

## Featured Trial Size Antibodies (10 µg)

Product Name	Cat. #	Species Reactivity	Applications
5-mC Monoclonal Antibody [33D3]	A-1014	Broad Range	DB, ELISA, IF, IHC, IP, MeDIP
5-hmC Monoclonal Antibody [HMC/4D9]	A-1018	Broad Range	DB, ELISA, IF, IHC, MeDIP
CRISPR/Cas9 Monoclonal Antibody [7A9]	A-9000	Species Independent	ELISA, IF, IP, WB
CRISPR/Cas9 (SaCas9) Monoclonal Antibody [6H4]	A-9001	Species Independent	ELISA, IF, IP, WB

## Trial Size Antibodies (20 µg)

Product Name	Cat. #	Species Reactivity	Applications
A4GNT Polyclonal Antibody	A62026	Human	ELISA, IHC
aac Polyclonal Antibody	A64338	Actinoplanes utahensis	ELISA
AADAC Polyclonal Antibody	A57918	Human	ELISA, IHC
AADACL2 Polyclonal Antibody	A62030	Human	ELISA, WB
AADAT Polyclonal Antibody	A66746	Human	ELISA, IHC
AANAT Polyclonal Antibody	A69907	Human	ELISA, WB, IF
AAO Polyclonal Antibody	A55231	Cucurbita maxima	ELISA, WB
AAP1 Polyclonal Antibody	A64314	Saccharomyces cerevisiae	ELISA
AAR2 Polyclonal Antibody	A57922	Human	ELISA, IHC
AARS Polyclonal Antibody	A68447	Human	ELISA, IHC, IF
AARSD1 Polyclonal Antibody	A57926	Human	ELISA, WB, IHC
AASDHPPT Polyclonal Antibody	A56783	Human	ELISA, WB, IHC
AATF Polyclonal Antibody	A51673	Human	ELISA, WB
ABCA12 Polyclonal Antibody	A69003	Human	ELISA, IF
ABCA9 Polyclonal Antibody	A66090	Human	ELISA, IF
ABCB1 Polyclonal Antibody	A52229	Human	ELISA, WB, IHC, IF
Abcb4 Polyclonal Antibody	A52844	Mouse	ELISA
ABCB5 Polyclonal Antibody	A57930	Human	ELISA, IHC, IF
ABCF3 Polyclonal Antibody	A57934	Human	ELISA, IHC
Abhd11 Polyclonal Antibody	A57938	Mouse	ELISA
ABHD14A Polyclonal Antibody	A57942	Human	ELISA, IHC, IF
ABHD14B Polyclonal Antibody	A57377	Human, Mouse	ELISA, WB, IHC
ABHD18 Polyclonal Antibody	A64610	Human	ELISA, IHC
ABHD2 Polyclonal Antibody	A57946	Human, Mouse, Rat	ELISA, WB, IHC, IF
ABHD3 Polyclonal Antibody	A51129	Human	ELISA, IHC, IF
ABI3 Polyclonal Antibody	A57950	Human, Mouse	ELISA, WB, IHC
ABI3BP Polyclonal Antibody	A65862	Human	ELISA, IHC
ABI5 Polyclonal Antibody	A54576	Arabidopsis thaliana	ELISA

Product Name	Cat. #	Species Reactivity	Applications
ABL1 Polyclonal Antibody	A51765	Human	ELISA
ABL2 Polyclonal Antibody	A69383	Human	ELISA, WB
ABLIM2 Polyclonal Antibody	A65550	Human	ELISA, IHC
ABLIM3 Polyclonal Antibody	A57954	Human	ELISA, IHC
ABTB1 Polyclonal Antibody	A57958	Human	ELISA
AC1 Polyclonal Antibody	A50768	African cassava mosaic virus	ELISA
ACAA1 Polyclonal Antibody	A50005	Human	ELISA, IHC
ACAD8 Polyclonal Antibody	A68138	Human, Mouse	ELISA, WB, IHC
ACADL Polyclonal Antibody	A57962	Human, Mouse	ELISA, WB, IHC
ACADM Polyclonal Antibody	A62038	Human	ELISA, IHC
ACAP1 Polyclonal Antibody	A64822	Human, Mouse	ELISA, WB, IHC
ACBD4 Polyclonal Antibody	A51045	Human, Mouse	ELISA, WB, IHC
ACBD5 Polyclonal Antibody	A50977	Human	ELISA, WB
ACBD7 Polyclonal Antibody	A57966	Human	ELISA
ACD Polyclonal Antibody	A70067	Human, Mouse	ELISA, WB
ACHE Polyclonal Antibody	A55190	Bovine	ELISA
ACKR1 Polyclonal Antibody	A50945	Human, Mouse	ELISA, WB, IF
ACLY Polyclonal Antibody	A62042	Human	ELISA, IHC, IF
ACOT11 Polyclonal Antibody	A57970	Human	ELISA
ACOT6 Polyclonal Antibody	A64966	Human	ELISA, IHC, IF
ACOT9 Polyclonal Antibody	A51433	Human	ELISA, IHC
ACOX3 Polyclonal Antibody	A62046	Human	ELISA
acpP Polyclonal Antibody	A56931	Staphylococcus aureus	ELISA
ACPP Polyclonal Antibody	A53618	Human	ELISA, IHC
acpS Polyclonal Antibody	A65374	Staphylococcus aureus	ELISA
ACRV1 Polyclonal Antibody	A57974	Human	ELISA, IHC
ACSBG1 Polyclonal Antibody	A67014	Human	ELISA, IHC
ACSBG2 Polyclonal Antibody	A57978	Human	ELISA, WB, IHC
ACSF2 Polyclonal Antibody	A66998	Human	ELISA, IHC
ACSL4 Polyclonal Antibody	A50628	Human	ELISA, WB, IHC
ACSL5 Polyclonal Antibody	A68074	Human, Rat	ELISA, WB, IHC
ACSL5 Polyclonal Antibody	A68078	Human	ELISA, IHC
Acta1 Polyclonal Antibody	A51958	Mouse, Rat, Zebrafish	ELISA, WB
ACTA1 Polyclonal Antibody	A54420	Human, Mouse, Rat	ELISA, WB, IHC
ACTA2 Polyclonal Antibody	A64156	Human	ELISA, IHC
ACTB Polyclonal Antibody	A52074	Human, Mouse	ELISA, WB, IHC, IF
Actb Polyclonal Antibody	A53186	Mouse, Human, Rat, Zebrafish	ELISA, WB
ACTG2 Polyclonal Antibody	A64164	Human	ELISA, WB, IHC
Actin, Alpha Skeletal Muscle Polyclonal Antibody	A57762	Human	ELISA
Actin, Alpha Skeletal Muscle Polyclonal Antibody	A57763	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
Actin, Cytoplasmic 1 Polyclonal Antibody	A57834	Human	ELISA, IHC, IF
ACTL7A Polyclonal Antibody	A54324	Human, Mouse	ELISA, WB, IHC
ACTL7B Polyclonal Antibody	A68234	Human, Mouse	ELISA, WB, IHC
ACTL8 Polyclonal Antibody	A57982	Human	ELISA, WB, IHC
ACTR10 Polyclonal Antibody	A57986	Human	ELISA, WB
ACTR8 Polyclonal Antibody	A57990	Human	ELISA, WB
ACTRT1 Polyclonal Antibody	A57994	Human	ELISA, IHC
ACTRT2 Polyclonal Antibody	A57998	Human	ELISA, WB
ACTRT3 Polyclonal Antibody	A58002	Human	ELISA, WB, IHC
ACVR1B Polyclonal Antibody	A58006	Human	ELISA
ADA Polyclonal Antibody	A50009	Human	ELISA, WB, IHC
ADAL Polyclonal Antibody	A65254	Human, Mouse	ELISA, WB
ADAM10 Polyclonal Antibody	A58010	Human	ELISA, IHC
ADAM15 Polyclonal Antibody	A52969	Human	ELISA
ADAM7 Polyclonal Antibody	A67258	Human	ELISA, IHC
ADAM8 Polyclonal Antibody	A69791	Human, Mouse	ELISA, WB, IHC, IF
ADAMDEC1 Polyclonal Antibody	A62050	Human	ELISA
ADAMTS12 Polyclonal Antibody	A62054	Human	ELISA, IHC
ADAMTS13 Polyclonal Antibody	A69387	Human	ELISA, IHC
ADAMTS15 Polyclonal Antibody	A70063	Human	ELISA, IHC
ADAMTS4 Polyclonal Antibody	A50013	Human, Mouse	ELISA, WB, IHC
ADAMTS7 Polyclonal Antibody	A58014	Human, Mouse	ELISA, WB, IHC
ADAP2 Polyclonal Antibody	A58018	Human	ELISA
ADAT1 Polyclonal Antibody	A58022	Human, Mouse	ELISA, WB, IHC
Adck2 Polyclonal Antibody	A54788	Mouse, Human	ELISA, WB, IHC
ADGRA3 Polyclonal Antibody	A66026	Human	ELISA, IF
ADGRL4 Polyclonal Antibody	A58026	Human	ELISA, WB, IHC, IF
ADH1B Polyclonal Antibody	A54784	Human	ELISA, WB, IHC
Adh5 Polyclonal Antibody	A55117	Mouse	ELISA, WB
ADH7 Polyclonal Antibody	A55635	Human, Mouse	ELISA, WB, IHC
Adh7 Polyclonal Antibody	A65186	Mouse	ELISA, WB
ADIPOQ Polyclonal Antibody	A53605	Human	ELISA, WB, IHC
ADIPOQ Polyclonal Antibody	A54069	Human, Mouse	ELISA, WB, IHC
ADIPOQ Polyclonal Antibody	A54866	bovine	ELISA
ADIPOQ Polyclonal Antibody	A55046	Bovine	ELISA, WB
ADIPOQ Polyclonal Antibody	A55747	Human	ELISA, IHC
Adipoq Polyclonal Antibody	A56646	Mouse	ELISA
ADM Polyclonal Antibody	A55846	Human	ELISA, IHC
ADO Polyclonal Antibody	A70075	Human	ELISA, IHC
ADORA1 Polyclonal Antibody	A62058	Human	ELISA, IF

Product Name	Cat. #	Species Reactivity	Applications
ADPRHL1 Polyclonal Antibody	A58030	Human	ELISA
Adrb2 Polyclonal Antibody	A57241	Rat	ELISA
ADRBK1 Polyclonal Antibody	A53488	Human	ELISA, IHC
AEN Polyclonal Antibody	A51057	Human	ELISA
AFG3L2 Polyclonal Antibody	A51605	Human	ELISA, WB, IHC
AFP Polyclonal Antibody	A52792	Human	ELISA, IHC
Afp Polyclonal Antibody	A55295	Mouse, Human, Rat	ELISA, WB
AFP Polyclonal Antibody	A55703	Human	ELISA
AFTPH Polyclonal Antibody	A62062	Human	ELISA, IHC
AGA Polyclonal Antibody	A56018	Human, Mouse	ELISA, WB
AGA2 Polyclonal Antibody	A50792	Saccharomyces cerevisiae	ELISA
AGBL2 Polyclonal Antibody	A58034	Human, Mouse	ELISA, WB, IHC
AGER Polyclonal Antibody	A52166	Human, Mouse	ELISA, IHC, IF
aglB Polyclonal Antibody	A54432	Aspergillus niger	ELISA
aglB Polyclonal Antibody	A55042	Aspergillus niger	ELISA
AGO2 Polyclonal Antibody	A68299	Human, Mouse	ELISA, WB, IHC, ChIP
AGPS Polyclonal Antibody	A69391	Human	ELISA, WB, IHC, IF
AGR3 Polyclonal Antibody	A58038	Human	ELISA
AGRN Polyclonal Antibody	A68451	Human	ELISA, IF
AGT Polyclonal Antibody	A50656	Human	ELISA, IHC, IF
Agxt Polyclonal Antibody	A50017	Mouse	ELISA
AGXT2 Polyclonal Antibody	A69195	Human	ELISA, WB, IF
AHCY Polyclonal Antibody	A54380	Human	ELISA, IHC
AHCY Polyclonal Antibody	A54894	Human	ELISA, IHC
AHNAK Polyclonal Antibody	A68819	Human	ELISA, IF
AHNAK2 Polyclonal Antibody	A70007	Human	ELISA, IHC, IF
AHSA1 Polyclonal Antibody	A51785	Human	ELISA, WB, IHC
AHSA2 Polyclonal Antibody	A65406	Human, Mouse, Rat	ELISA, WB, IHC
AHSG Polyclonal Antibody	A52102	Human	ELISA, IHC, IF
AIF1 Polyclonal Antibody	A55283	Human	ELISA, IHC, IF
Aif1 Polyclonal Antibody	A56194	Mouse, Human	ELISA, WB
Aif1 Polyclonal Antibody	A62066	Rat	ELISA
AIFM1 Polyclonal Antibody	A56759	Human	ELISA
aiiM Polyclonal Antibody	A64582	Microbacterium testaceum	ELISA
AIMP1 Polyclonal Antibody	A69867	Human	ELISA, WB, IHC, IF
AK1 Polyclonal Antibody	A50021	Human	ELISA, IHC
AKAP7 Polyclonal Antibody	A58042	Human	ELISA, IHC
AKR1C2 Polyclonal Antibody	A56519	Human, Rat	ELISA, WB
AKR1D1 Polyclonal Antibody	A62070	Human	ELISA
AKT1 Polyclonal Antibody	A68300	Human	ELISA, WB, IHC, ChIP

Product Name	Cat. #	Species Reactivity	Applications
AKT2 Polyclonal Antibody	A56166	Human	ELISA, IHC
AKTIP Polyclonal Antibody	A57353	Human	ELISA
alaC Polyclonal Antibody	A55431	Escherichia coli	ELISA
ALAS2 Polyclonal Antibody	A62074	Human	ELISA, WB, IHC
ALB Polyclonal Antibody	A55898	Dog	ELISA
ALB Polyclonal Antibody	A56619	Macaca fascicularis	ELISA
Aldh1a1 Polyclonal Antibody	A58046	Mouse	ELISA, WB
ALDH1L2 Polyclonal Antibody	A64914	Human	ELISA, IHC
ALDH3A1 Polyclonal Antibody	A50025	Human	ELISA, WB, IHC
ALDH3B2 Polyclonal Antibody	A50029	Human	ELISA, IHC
ALDH5A1 Polyclonal Antibody	A57337	Human, Mouse, Rat	ELISA, WB, IHC
ALDH7A1 Polyclonal Antibody	A68455	Human, Mouse, Rat	ELISA, WB, IF
ALDOA Polyclonal Antibody	A52360	Human, Mouse	ELISA, WB, IHC
ALG12 Polyclonal Antibody	A67914	Human	ELISA, IHC, IF
ALG14 Polyclonal Antibody	A58050	Human	ELISA, IHC
ALG5 Polyclonal Antibody	A58054	Human	ELISA, IHC
ALG9 Polyclonal Antibody	A69203	Human, Rat	ELISA, WB
ALKBH3 Polyclonal Antibody	A57365	Human	ELISA, IHC, IF
Allergen Api g 1.0201 Polyclonal Antibody	A50728	Apium graveolens	ELISA
Allergen Ara h 1 Polyclonal Antibody	A64330	Peanut	ELISA
Allergen Car b 1 Polyclonal Antibody	A57393	Carpinus betulus	ELISA, WB
Allergen Dac g 3 Polyclonal Antibody	A64242	Dactylis glomerata	ELISA
Allergen Dac g 4 Polyclonal Antibody	A64230	Dactylis glomerata	ELISA
ALMS1P Polyclonal Antibody	A58058	Human	ELISA
ALOX15B Polyclonal Antibody	A56791	Human	ELISA, WB
ALOX5 Polyclonal Antibody	A58062	Human	ELISA, IHC
Alpha-1-Acid Glycoprotein 1 Polyclonal Antibody	A57652	Human	ELISA
Alpha1-Acidglycoprotein Polyclonal Antibody	A53849	Human	ELISA, WB
Alpha-insect Toxin LqhaIT Polyclonal Antibody	A69815	Leiurus quinquestriatus hebraeus	ELISA
Alpha-TCS Polyclonal Antibody	A64470	Trichosanthes kirilowii	ELISA
ALPI Polyclonal Antibody	A52182	Human	ELISA, WB
ALPL Polyclonal Antibody	A53198	Mouse	ELISA
ALPL Polyclonal Antibody	A54165	Human	ELISA, IHC
ALS2CR11 Polyclonal Antibody	A65042	Human	ELISA, IHC
ALS2CR12 Polyclonal Antibody	A69139	Human	ELISA, WB, IF
ALTA1 Polyclonal Antibody	A57453	Alternaria alternata	ELISA
AMA-1 Polyclonal Antibody	A53015	Plasmodium falciparum	ELISA
Amatoxin Polyclonal Antibody	A64220	Amanita phalloides	ELISA
Amatoxin Polyclonal Antibody	A64221	Amanita phalloides	ELISA
AMBN Polyclonal Antibody	A70287	Human, Rat	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
AMBP Polyclonal Antibody	A52265	Human	ELISA, IF
AMBP Polyclonal Antibody	A53628	Human, Mouse	ELISA, WB, IHC
AMDHD1 Polyclonal Antibody	A55982	Human	ELISA
Amh Polyclonal Antibody	A54922	Mouse, Human	ELISA, WB, IHC
AMH Polyclonal Antibody	A56138	Human, Mouse	ELISA, WB, IHC, IF
amiF Polyclonal Antibody	A64558	Bradyrhizobium sp.	ELISA
AMIGO1 Polyclonal Antibody	A65886	Human	ELISA, IHC
AMMECR1 Polyclonal Antibody	A68771	Human	ELISA, IHC, IF
ampC Polyclonal Antibody	A56915	Pseudomonas aeruginosa	ELISA
AMTN Polyclonal Antibody	A62078	Human	ELISA, IHC
Amyloid Beta A4 Protein A Polyclonal Antibody	A57755	Human	ELISA, WB
Amyloid Beta A4 Protein A Polyclonal Antibody	A57756	Human, Mouse	ELISA, WB
ANAPC5 Polyclonal Antibody	A53997	Human	ELISA, IHC
ANG Polyclonal Antibody	A52942	Human	ELISA, IHC
ANG1 Polyclonal Antibody	A57654	Bovine	ELISA
ANGPT1 Polyclonal Antibody	A57770	Human	ELISA, WB, IHC
ANGPT2 Polyclonal Antibody	A52448	Human	ELISA, IHC
Angpt2 Polyclonal Antibody	A52516	Mouse	ELISA, WB, IHC
ANGPTL2 Polyclonal Antibody	A51389	Human, Mouse	ELISA, WB, IHC
ANGPTL2 Polyclonal Antibody	A51393	Human	ELISA, IHC
ANGPTL3 Polyclonal Antibody	A64080	Human	ELISA, IHC
ANGPTL7 Polyclonal Antibody	A50033	Human	ELISA, WB, IHC
ANGPTL8 Polyclonal Antibody	A54736	Human, Mouse	ELISA, WB, IHC
Angptl8 Polyclonal Antibody	A55287	Mouse	ELISA
ANK3 Polyclonal Antibody	A69875	Human	ELISA, IHC, IF
ANKDD1A Polyclonal Antibody	A64982	Human	ELISA, IHC
ANKMY1 Polyclonal Antibody	A58066	Human	ELISA, IHC
ANKMY2 Polyclonal Antibody	A58070	Human	ELISA, IHC
ANKRA2 Polyclonal Antibody	A67510	Human	ELISA, IHC
ANKRD10 Polyclonal Antibody	A58074	Human, Mouse	ELISA, WB
ANKRD23 Polyclonal Antibody	A65982	Human, Mouse	ELISA, WB, IHC
ANKRD33 Polyclonal Antibody	A65790	Human	ELISA, IHC
ANKRD49 Polyclonal Antibody	A66554	Human	ELISA, WB, IHC
ANKRD55 Polyclonal Antibody	A64918	Human	ELISA, IHC, IF
ANKRD7 Polyclonal Antibody	A58078	Human	ELISA
Annexin A5 Polyclonal Antibody	A50037	Cynops pyrrhogaster	ELISA
ANO2 Polyclonal Antibody	A69283	Human	ELISA, IHC, IF
ANO4 Polyclonal Antibody	A64886	Human	ELISA, IHC, IF
ANXA1 Polyclonal Antibody	A52396	Human	ELISA, WB
Anxa1 Polyclonal Antibody	A53424	Mouse, Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
Anxa1 Polyclonal Antibody	A53550	Mouse, Human	ELISA, WB
ANXA2 Polyclonal Antibody	A52816	Human	ELISA, WB
ANXA2 Polyclonal Antibody	A56863	Human, Mouse	ELISA, WB, IHC, IF, IP
ANXA3 Polyclonal Antibody	A52352	Human	ELISA
ANXA5 Polyclonal Antibody	A51897	Human, Mouse	ELISA, WB, IHC
ANXA5 Polyclonal Antibody	A56414	Human	ELISA, WB, IHC
ANXA6 Polyclonal Antibody	A53041	Human, Rat	ELISA, WB
ANXA9 Polyclonal Antibody	A55711	Human, Mouse	ELISA, WB, IHC
AOX1 Polyclonal Antibody	A53444	Human	ELISA, IHC
AP1S1 Polyclonal Antibody	A58082	Human	ELISA
AP1S3 Polyclonal Antibody	A53257	Human	ELISA, WB
AP24 Polyclonal Antibody	A64274	Nicotiana tabacum	ELISA, WB
AP2M1 Polyclonal Antibody	A53865	Human	ELISA, WB, IHC
AP2M1 Polyclonal Antibody	A54500	Human, Mouse, Rat	ELISA, WB, IHC
AP2S1 Polyclonal Antibody	A51809	Human	ELISA, IHC, IF
AP3M2 Polyclonal Antibody	A62082	Human	ELISA, IHC
AP3S2 Polyclonal Antibody	A58086	Human	ELISA, WB
AP4B1 Polyclonal Antibody	A58090	Human	ELISA, WB, IHC
APAF1 Polyclonal Antibody	A58094	Human	ELISA, IHC
apaLIM Polyclonal Antibody	A50869	Acetobacter pasteurianus	ELISA
APEH Polyclonal Antibody	A62086	Human, Mouse	ELISA, WB, IHC, IF
Apelin Polyclonal Antibody	A57832	Human	ELISA, IHC
Apelin Polyclonal Antibody	A57833	Human	ELISA, IHC
APEX1 Polyclonal Antibody	A68301	Human, Mouse	ELISA, WB, IHC, ChIP
APLNR Polyclonal Antibody	A58098	Human	ELISA, IHC, IF
APOA1 Polyclonal Antibody	A54336	Human	ELISA
APOA1 Polyclonal Antibody	A55259	Bovine, Mouse	ELISA, WB
Apoa4 Polyclonal Antibody	A53040	Mouse	ELISA
APOA4 Polyclonal Antibody	A57173	Human	ELISA, WB
APOB Polyclonal Antibody	A58102	Human	ELISA
APOBEC1 Polyclonal Antibody	A69395	Human	ELISA, IF
APOBEC3A Polyclonal Antibody	A58106	Human	ELISA
APOC3 Polyclonal Antibody	A55307	Human, Mouse, Rat	ELISA, WB, IHC
APOD Polyclonal Antibody	A50041	Human	ELISA, IHC
Apoe Polyclonal Antibody	A53280	Mouse	ELISA, WB
APOOL Polyclonal Antibody	A62090	Human	ELISA, IHC
Apoptin Polyclonal Antibody	A57857	Chicken anemia virus	ELISA, WB
APP Polyclonal Antibody	A54049	Human	ELISA, IHC
APP Polyclonal Antibody	A56146	Human, Mouse	ELISA, WB, IHC, IF
APPL1 Polyclonal Antibody	A68030	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
APX1 Polyclonal Antibody	A52772	Arabidopsis thaliana	ELISA, WB
APX2 Polyclonal Antibody	A52937	Arabidopsis thaliana	ELISA, WB
AQP1 Polyclonal Antibody	A53228	Human, Mouse, Rat	ELISA, WB, IF
AQP10 Polyclonal Antibody	A69135	Human	ELISA, IF
AQP4 Polyclonal Antibody	A50672	Human, Mouse	ELISA, WB
Arachin Ahy-3 Polyclonal Antibody	A56270	Peanut	ELISA
ARF4 Polyclonal Antibody	A52593	Human	ELISA, WB, IHC
ARF5 Polyclonal Antibody	A51477	Human	ELISA
ARF6 Polyclonal Antibody	A56150	Human, Rat	ELISA, WB, IHC
ARFIP2 Polyclonal Antibody	A58110	Human, Mouse	ELISA, WB, IHC
ARFRP1 Polyclonal Antibody	A69891	Human	ELISA, IHC
Arg1 Polyclonal Antibody	A58114	Mouse	ELISA
Arg2 Polyclonal Antibody	A58118	Mouse, Human	ELISA, WB
argF Polyclonal Antibody	A54173	Escherichia coli (strain K12)	ELISA
ARHGAP11A Polyclonal Antibody	A65518	Human	ELISA
ARHGAP19 Polyclonal Antibody	A55427	Human	ELISA, IHC
ARHGAP25 Polyclonal Antibody	A58122	Human	ELISA, IHC
ARHGAP27 Polyclonal Antibody	A65402	Human	ELISA, IHC
ARHGAP29 Polyclonal Antibody	A58126	Human	ELISA
ARHGDI A Polyclonal Antibody	A52636	Human	ELISA, WB, IHC
ARHGFE18 Polyclonal Antibody	A51713	Human	ELISA, WB
ARHGFE18 Polyclonal Antibody	A63334	Human	ELISA
ARHGFE39 Polyclonal Antibody	A58130	Human	ELISA
ARHGFE7 Polyclonal Antibody	A53117	Human	ELISA, WB
ARL1 Polyclonal Antibody	A68459	Human	ELISA, IHC, IF
ARL11 Polyclonal Antibody	A58134	Human	ELISA, IHC
ARL14 Polyclonal Antibody	A58138	Human	ELISA, WB, IHC
ARL15 Polyclonal Antibody	A67754	Human	ELISA, IHC
ARL16 Polyclonal Antibody	A64626	Human	ELISA, IHC
ARL2 Polyclonal Antibody	A52648	Human	ELISA, WB
ARL4A Polyclonal Antibody	A50045	Human	ELISA
ARL4C Polyclonal Antibody	A62094	Human	ELISA, IHC
ARL4D Polyclonal Antibody	A57149	Human	ELISA, IHC
ARL6IP1 Polyclonal Antibody	A64686	Human	ELISA, IHC, IF
ARL6IP5 Polyclonal Antibody	A62098	Human	ELISA, IHC, IF
ARL8A Polyclonal Antibody	A66682	Human	ELISA
ARL8B Polyclonal Antibody	A58142	Human, Mouse	ELISA, WB, IHC
ARMC10 Polyclonal Antibody	A58146	Human, Mouse	ELISA, WB, IHC
ARMC7 Polyclonal Antibody	A58150	Human	ELISA, IHC
ARMC8 Polyclonal Antibody	A58154	Human	ELISA



Product Name	Cat. #	Species Reactivity	Applications
ARMC9 Polyclonal Antibody	A58158	Human	ELISA, WB, IHC
ARMCX1 Polyclonal Antibody	A58162	Human	ELISA, IHC
ARMCX2 Polyclonal Antibody	A58166	Human, Mouse	ELISA, WB, IHC
ARMCX6 Polyclonal Antibody	A58170	Human	ELISA, WB, IHC
ARMT1 Polyclonal Antibody	A67382	Human	ELISA, WB, IHC
aroD Polyclonal Antibody	A50784	Salmonella typhi	ELISA
aroD Polyclonal Antibody	A57449	Salmonella typhi	ELISA
ARPC2 Polyclonal Antibody	A51677	Human	ELISA
ARPC3 Polyclonal Antibody	A51709	Human	ELISA, WB, IHC, IF
ARPIN Polyclonal Antibody	A65910	Human	ELISA, WB, IHC
ARRB1 Polyclonal Antibody	A62102	Human	ELISA
ARRDC1-AS1 Polyclonal Antibody	A67946	Human	ELISA, IHC
ARRDC3 Polyclonal Antibody	A58174	Human, Mouse	ELISA, WB
ARSB Polyclonal Antibody	A54292	Human, Mouse	ELISA, WB
ARSE Polyclonal Antibody	A69399	Human	ELISA, IHC, IF
ARSG Polyclonal Antibody	A50049	Human	ELISA, WB, IHC
ARSH Polyclonal Antibody	A68915	Human	ELISA, WB, IF
ARSI Polyclonal Antibody	A65054	Human	ELISA, IHC
ART5 Polyclonal Antibody	A66466	Human	ELISA, WB, IHC
ARV1 Polyclonal Antibody	A51277	Human	ELISA, IF
AS3MT Polyclonal Antibody	A69299	Human	ELISA, IHC, IF
ASAH2 Polyclonal Antibody	A62106	Human, Rat	ELISA, WB, IHC
ASAP3 Polyclonal Antibody	A66542	Human	ELISA, WB, IHC
ASB10 Polyclonal Antibody	A66798	Human	ELISA, IHC, IF
ASB11 Polyclonal Antibody	A67066	Human	ELISA, IHC
ASB13 Polyclonal Antibody	A58178	Human	ELISA, IHC
ASB6 Polyclonal Antibody	A58182	Human	ELISA, IHC
ASB7 Polyclonal Antibody	A67274	Human	ELISA, WB, IHC
ASB8 Polyclonal Antibody	A58186	Human	ELISA, WB, IHC
ASB9 Polyclonal Antibody	A54296	Human	ELISA
ASCC1 Polyclonal Antibody	A58190	Human, Mouse	ELISA, WB, IHC
ASF1B Polyclonal Antibody	A67618	Human	ELISA, WB, IHC
ASGR1 Polyclonal Antibody	A52388	Human	ELISA
ASGR2 Polyclonal Antibody	A62110	Human	ELISA, IHC, IF
ASH2L Polyclonal Antibody	A56302	Human, Mouse	ELISA, WB, IHC, IF
ASIC2 Polyclonal Antibody	A69851	Human	ELISA, IHC, IF
ASL Polyclonal Antibody	A62114	Human	ELISA, IHC
ASMT Polyclonal Antibody	A68463	Human	ELISA, IF
ASNSD1 Polyclonal Antibody	A67866	Human	ELISA, IHC
ASPH Polyclonal Antibody	A54361	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
ASPHD1 Polyclonal Antibody	A65182	Human	ELISA, IHC
ASPHD2 Polyclonal Antibody	A58194	Human	ELISA, WB
ASTL Polyclonal Antibody	A65638	Human	ELISA, IHC
ASZ1 Polyclonal Antibody	A68286	Human, Rat	ELISA, WB, IHC, IF
At2g24580 Polyclonal Antibody	A53392	Arabidopsis thaliana	ELISA, WB
ATAD3C Polyclonal Antibody	A65174	Human	ELISA, IHC
ATAT1 Polyclonal Antibody	A68935	Human, Mouse, Rat	ELISA, WB, IHC
ATF3 Polyclonal Antibody	A62118	Human	ELISA, IHC
ATG1 Polyclonal Antibody	A65366	Candida glabrata	ELISA
ATG10 Polyclonal Antibody	A67182	Human	ELISA, IF
ATG16L1 Polyclonal Antibody	A50989	Human	ELISA, IHC
ATG2B Polyclonal Antibody	A58198	Human	ELISA, IHC
ATG3 Polyclonal Antibody	A55479	Human	ELISA, WB, IHC
ATG4A Polyclonal Antibody	A51133	Human, Mouse	ELISA, WB
ATG4D Polyclonal Antibody	A65918	Human, Mouse	ELISA, WB, IHC, IF
ATG5 Polyclonal Antibody	A64172	Human	ELISA, IHC, IF
ATG7 Polyclonal Antibody	A62122	Human	ELISA, WB, IHC, IF
ATG8 Polyclonal Antibody	A65370	Candida glabrata	ELISA
ATIC Polyclonal Antibody	A62126	Human	ELISA, IHC
ATL2 Polyclonal Antibody	A66526	Human	ELISA, WB, IHC, IF
ATL3 Polyclonal Antibody	A50981	Human	ELISA, WB, IHC, IF
ATOH1 Polyclonal Antibody	A56014	Human	ELISA, IHC
Atoh1 Polyclonal Antibody	A62130	Mouse	ELISA, WB
ATP12A Polyclonal Antibody	A62134	Human	ELISA, IHC, IF
ATP1B3 Polyclonal Antibody	A58202	Human	ELISA, IHC
ATP2A2 Polyclonal Antibody	A58206	Human, Mouse	ELISA, WB, IHC, IF
ATP4B Polyclonal Antibody	A53516	Human	ELISA
ATP5A1 Polyclonal Antibody	A50061	Human	ELISA, WB
ATP5A1 Polyclonal Antibody	A52660	Human	ELISA, WB, IHC
Atp5b Polyclonal Antibody	A54966	Mouse	ELISA
ATP5D Polyclonal Antibody	A50065	Human	ELISA, IHC
ATP5I Polyclonal Antibody	A58210	Human	ELISA
ATP6AP2 Polyclonal Antibody	A57277	Human, Mouse	ELISA, WB
ATP6V0A2 Polyclonal Antibody	A68222	Human	ELISA, IF
ATP6V1E1 Polyclonal Antibody	A52724	Human	ELISA, WB, IF
ATP6V1E2 Polyclonal Antibody	A58214	Human	ELISA
ATP6V1F Polyclonal Antibody	A52952	Human	ELISA, IHC, IF
ATP6V1G1 Polyclonal Antibody	A51613	Human	ELISA, WB, IHC
ATP6V1G2 Polyclonal Antibody	A62138	Human, Mouse, Rat	ELISA, WB, IHC
ATPAF1 Polyclonal Antibody	A69403	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
ATPAF2 Polyclonal Antibody	A54269	Human	ELISA
atpB Polyclonal Antibody	A53460	Lemna minor	ELISA
atpE Polyclonal Antibody	A57061	Strain OhILAs	ELISA
ATXN3 Polyclonal Antibody	A50069	Human, Mouse	ELISA, WB, IHC
ATXN7L1 Polyclonal Antibody	A68094	Human	ELISA, IHC
ATXN7L3 Polyclonal Antibody	A58218	Human	ELISA, IHC
AURKB Polyclonal Antibody	A50073	Human	ELISA, IHC
AVP Polyclonal Antibody	A54862	Human, Rat	ELISA, WB, IHC
AVPI1 Polyclonal Antibody	A65322	Human	ELISA, IHC
Avpr1a Polyclonal Antibody	A53272	Rat, Mouse, Human	ELISA, WB, IHC
Avpr1b Polyclonal Antibody	A53413	Rat	ELISA, WB
AZGP1 Polyclonal Antibody	A58222	Human, Mouse	ELISA, WB, IHC, IP
AZIN1 Polyclonal Antibody	A62142	Human, Mouse	ELISA, WB, IHC
B2M Polyclonal Antibody	A57870	Human, Mouse	ELISA, WB, IHC
B3GALT2 Polyclonal Antibody	A58226	Human	ELISA, IHC
B3GNT2 Polyclonal Antibody	A58230	Human	ELISA, WB, IHC
B3GNT5 Polyclonal Antibody	A50081	Human	ELISA, IHC
B4GALNT2 Polyclonal Antibody	A67050	Human	ELISA, IHC
B4GALT3 Polyclonal Antibody	A53869	Human	ELISA, WB, IHC
B4GALT5 Polyclonal Antibody	A64594	Human, Mouse	ELISA, WB, IHC
B4GAT1 Polyclonal Antibody	A50077	Human	ELISA, IHC
BABAM2 Polyclonal Antibody	A70247	Human	ELISA, IHC, IF
Bag3 Polyclonal Antibody	A55866	Mouse, Human	ELISA, WB
BAG5 Polyclonal Antibody	A55319	Human, Mouse	ELISA, WB, IHC
BAGE3 Polyclonal Antibody	A65774	Human	ELISA, IHC
BAIAP2L1 Polyclonal Antibody	A58234	Human	ELISA, IHC
BAIAP3 Polyclonal Antibody	A68467	Human	ELISA, IHC
BALF5 Polyclonal Antibody	A56903	Epstein-Barr virus	ELISA
bar Polyclonal Antibody	A54543	Streptomyces hygroscopicus	ELISA
BARX1 Polyclonal Antibody	A58238	Human	ELISA, WB, IHC
BATF Polyclonal Antibody	A58242	Human, Mouse	ELISA, WB, IHC
BATF2 Polyclonal Antibody	A66342	Human	ELISA, IHC
Batf3 Polyclonal Antibody	A50085	Rat	ELISA
BAZ2B Polyclonal Antibody	A69271	Human	ELISA, IF
BBE1 Polyclonal Antibody	A54365	Eschscholzia californica	ELISA
BBS7 Polyclonal Antibody	A70011	Human	ELISA, IF
BCAP31 Polyclonal Antibody	A51753	Human	ELISA, WB, IHC, IF
BCAT1 Polyclonal Antibody	A62146	Human	ELISA, IHC, IF
BCAT2 Polyclonal Antibody	A55902	Human, Mouse	ELISA, WB, IHC
BCDIN3D Polyclonal Antibody	A65746	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
BCHE Polyclonal Antibody	A53245	Dog	ELISA, IHC
BCKDHA Polyclonal Antibody	A69407	Human	ELISA, WB, IHC
BCL10 Polyclonal Antibody	A55074	Human	ELISA, IHC
Bcl2 Polyclonal Antibody	A52548	Mouse, Human	ELISA, IHC
BCL2L1 Polyclonal Antibody	A56459	Human	ELISA
BCL2L2 Polyclonal Antibody	A56350	Human	ELISA, IHC
BCO1 Polyclonal Antibody	A67286	Human	ELISA, IHC
BCOR Polyclonal Antibody	A65658	Human	ELISA
BDH1 Polyclonal Antibody	A53205	Human, Mouse	ELISA, WB, IHC
BDH2 Polyclonal Antibody	A58246	Human	ELISA, IHC
BDNF Polyclonal Antibody	A52601	Human, Mouse	ELISA, WB, IHC
BDNF Polyclonal Antibody	A53065	Rat	ELISA, WB
BDNF Polyclonal Antibody	A58250	Human	ELISA, IHC
BEND7 Polyclonal Antibody	A58254	Human	ELISA, IHC
Beta-mammal Toxin Css4 Polyclonal Antibody	A69827	Centruroides suffusus	ELISA
BEX1 Polyclonal Antibody	A67210	Human	ELISA, IHC
BEX2 Polyclonal Antibody	A58258	Human	ELISA, IHC
BEX5 Polyclonal Antibody	A65114	Human	ELISA, IHC
BGLAP Polyclonal Antibody	A52249	Human	ELISA, IHC
Bglap Polyclonal Antibody	A57033	Rat	ELISA
BGN Polyclonal Antibody	A62150	Human	ELISA, IHC, IF
BHLHE41 Polyclonal Antibody	A69275	Human	ELISA, WB
BHMT2 Polyclonal Antibody	A64140	Human, Mouse	ELISA, WB, IHC, IF
BID Polyclonal Antibody	A55679	Human	ELISA
BIN2 Polyclonal Antibody	A68471	Human	ELISA, IF
Birc5 Polyclonal Antibody	A56487	Mouse	ELISA
bla Polyclonal Antibody	A56378	Escherichia coli	ELISA
blaNDM-1 Polyclonal Antibody	A55774	Klebsiella pneumoniae	ELISA
blaNDM-1 Polyclonal Antibody	A64578	Klebsiella pneumoniae	ELISA
BLLF1 Polyclonal Antibody	A64514	Epstein-Barr virus	ELISA
BLOC1S1 Polyclonal Antibody	A53861	Human	ELISA, WB, IHC, IF
BLOC1S5 Polyclonal Antibody	A66538	Human	ELISA, IHC
BMI1 Polyclonal Antibody	A53945	Human	ELISA, WB, IHC
BMP2 Polyclonal Antibody	A54951	Human	ELISA, WB, IHC, IF
BMP3 Polyclonal Antibody	A52753	Human, Mouse	ELISA, WB, IHC
BMP4 Polyclonal Antibody	A55459	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC
BMP6 Polyclonal Antibody	A52078	Human	ELISA, IHC
BMP7 Polyclonal Antibody	A52444	Human, Mouse	ELISA, WB, IHC, IF
BMPER Polyclonal Antibody	A70079	Human	ELISA, IHC
BMPR1A Polyclonal Antibody	A50089	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
BOLA1 Polyclonal Antibody	A62154	Human	ELISA
botA Polyclonal Antibody	A57457	Clostridium botulinum	ELISA
botB Polyclonal Antibody	A50736	Clostridium botulinum	ELISA
botD Polyclonal Antibody	A50764	Clostridium botulinum	ELISA
Botulinum Neurotoxin Type C1 Polyclonal Antibody	A50780	Clostridium botulinum	ELISA
Botulinum Neurotoxin Type E Polyclonal Antibody	A57517	Clostridium butyricum	ELISA
Bovine Serum Albumin Polyclonal Antibody	A55766	Bovine	ELISA
BRAP Polyclonal Antibody	A69979	Human, Rat	ELISA, WB, IHC, IF
BRCA1 Polyclonal Antibody	A54101	Human	ELISA
BRCC3 Polyclonal Antibody	A54460	Human	ELISA, WB, IHC
BRINP1 Polyclonal Antibody	A62382	Human	ELISA, IHC, IF
BRINP2 Polyclonal Antibody	A58262	Human	ELISA, IHC
BSG Polyclonal Antibody	A52086	Human	ELISA, WB, IHC
BTBD2 Polyclonal Antibody	A53993	Human	ELISA, WB, IHC, IF
BTD Polyclonal Antibody	A69411	Human	ELISA, IHC
BTF3 Polyclonal Antibody	A51857	Human	ELISA, IHC, IF
BTG4 Polyclonal Antibody	A67634	Human, Mouse	ELISA, WB, IHC
BTN2A1 Polyclonal Antibody	A58266	Human	ELISA, IHC
BTN2A2 Polyclonal Antibody	A51149	Human, Mouse	ELISA, WB, IHC
BTN3A1 Polyclonal Antibody	A58270	Human, Mouse	ELISA, WB, IHC
BTN3A2 Polyclonal Antibody	A58274	Human	ELISA
BTN3A3 Polyclonal Antibody	A58278	Human	ELISA
BTNL3 Polyclonal Antibody	A65718	Human	ELISA, IHC
BTNL9 Polyclonal Antibody	A65438	Human	ELISA
BUD31 Polyclonal Antibody	A63098	Human, Mouse	ELISA, WB, IHC, IF
BZLF2 Polyclonal Antibody	A69831	Epstein-Barr virus	ELISA, WB
C10orf111 Polyclonal Antibody	A58282	Human	ELISA, IHC, IF
C10orf32 Polyclonal Antibody	A66678	Human	ELISA, IHC, IF
C10orf53 Polyclonal Antibody	A58286	Human	ELISA, IHC
C10orf54 Polyclonal Antibody	A67834	Human, Mouse	ELISA, WB, IHC
C10orf62 Polyclonal Antibody	A58290	Human	ELISA, IHC
C10orf82 Polyclonal Antibody	A66558	Human, Mouse	ELISA, WB, IHC
C10orf91 Polyclonal Antibody	A65230	Human	ELISA, IHC
C11orf1 Polyclonal Antibody	A67706	Human	ELISA, IHC
C11orf45 Polyclonal Antibody	A58294	Human	ELISA, IHC
C11orf53 Polyclonal Antibody	A66038	Human, Mouse	ELISA, WB, IHC
C11orf68 Polyclonal Antibody	A58298	Human, Mouse	ELISA, WB, IHC
C11orf73 Polyclonal Antibody	A50969	Human	ELISA, IHC
C11orf84 Polyclonal Antibody	A67910	Human	ELISA
C11orf85 Polyclonal Antibody	A64902	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
C12orf43 Polyclonal Antibody	A66910	Human	ELISA, IHC
C12orf49 Polyclonal Antibody	A51281	Human	ELISA, IHC
C12orf50 Polyclonal Antibody	A66386	Human	ELISA
C12orf57 Polyclonal Antibody	A58302	Human	ELISA
C12orf60 Polyclonal Antibody	A58306	Human	ELISA, IHC, IF
C15orf32 Polyclonal Antibody	A64898	Human	ELISA
C15orf40 Polyclonal Antibody	A58310	Human	ELISA, IHC
C15orf41 Polyclonal Antibody	A58314	Human	ELISA, WB
C15orf53 Polyclonal Antibody	A66962	Human	ELISA
C15orf57 Polyclonal Antibody	A67790	Human, Mouse	ELISA, WB
C16orf78 Polyclonal Antibody	A58318	Human	ELISA, IHC
C16orf87 Polyclonal Antibody	A65702	Human	ELISA, WB
C17orf47 Polyclonal Antibody	A66970	Human	ELISA, IHC
C17orf58 Polyclonal Antibody	A64858	Human	ELISA, IHC, IF
C17orf62 Polyclonal Antibody	A67150	Human, Mouse	ELISA, WB, IHC, IF
C17orf64 Polyclonal Antibody	A65850	Human, Mouse	ELISA, WB, IHC
C17orf75 Polyclonal Antibody	A67718	Human	ELISA
C17orf78 Polyclonal Antibody	A66854	Human	ELISA, IHC, IF
C19orf44 Polyclonal Antibody	A58322	Human	ELISA, IHC
C19orf47 Polyclonal Antibody	A58326	Human, Mouse	ELISA, WB, IHC
C19orf53 Polyclonal Antibody	A58330	Human	ELISA
C1D Polyclonal Antibody	A56635	Human	ELISA, IHC
C1GALT1 Polyclonal Antibody	A70295	Human	ELISA, WB, IHC, IF
C1orf101 Polyclonal Antibody	A65202	Human, Mouse	ELISA, WB, IHC, IF
C1orf105 Polyclonal Antibody	A62170	Human	ELISA, IHC
C1orf111 Polyclonal Antibody	A58334	Human	ELISA, IHC
C1orf112 Polyclonal Antibody	A67606	Human	ELISA, IHC
C1orf123 Polyclonal Antibody	A58338	Human, Mouse	ELISA, WB, IHC
C1orf127 Polyclonal Antibody	A67046	Human	ELISA, IHC
C1orf158 Polyclonal Antibody	A58342	Human	ELISA
C1orf189 Polyclonal Antibody	A65286	Human	ELISA, IHC, IF
C1orf210 Polyclonal Antibody	A66022	Human	ELISA
C1orf53 Polyclonal Antibody	A65210	Human	ELISA, IHC
C1orf54 Polyclonal Antibody	A58346	Human	ELISA, IHC
C1orf64 Polyclonal Antibody	A66770	Human	ELISA
C1orf87 Polyclonal Antibody	A66838	Human	ELISA
C1QTNF1 Polyclonal Antibody	A51293	Human	ELISA, IHC
C1QTNF3 Polyclonal Antibody	A52873	Human	ELISA, IHC
C1QTNF6 Polyclonal Antibody	A67546	Human, Mouse	ELISA, WB, IHC
C20orf195 Polyclonal Antibody	A58350	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
C20orf196 Polyclonal Antibody	A66102	Human	ELISA
C20orf197 Polyclonal Antibody	A66346	Human	ELISA
C20orf27 Polyclonal Antibody	A58354	Human	ELISA, WB, IHC
C21orf2 Polyclonal Antibody	A62174	Human	ELISA
C21orf33 Polyclonal Antibody	A53901	Human	ELISA, WB, IHC
C21orf62-AS1 Polyclonal Antibody	A64662	Human	ELISA
C22orf23 Polyclonal Antibody	A67462	Human	ELISA
C2orf15 Polyclonal Antibody	A58358	Human	ELISA, WB, IHC
C2orf47 Polyclonal Antibody	A51197	Human, Mouse	ELISA, WB, IHC
C2orf49 Polyclonal Antibody	A67178	Human	ELISA, IHC
C2orf50 Polyclonal Antibody	A66610	Human	ELISA
C2orf69 Polyclonal Antibody	A58362	Human	ELISA, IHC
C2orf76 Polyclonal Antibody	A64906	Human	ELISA, IHC
C2orf80 Polyclonal Antibody	A64598	Human	ELISA
C2orf83 Polyclonal Antibody	A65142	Human	ELISA, IHC
C3 Polyclonal Antibody	A52432	Human	ELISA, IHC, IF
C3 Polyclonal Antibody	A55275	Mouse, Human	ELISA, WB
C3orf14 Polyclonal Antibody	A67846	Human, Mouse, Rat	ELISA, WB, IHC
C3orf18 Polyclonal Antibody	A68130	Human, Mouse	ELISA, WB, IHC, IF
C3orf20 Polyclonal Antibody	A58366	Human	ELISA, IF
C3orf36 Polyclonal Antibody	A64938	Human	ELISA, IHC
C3orf49 Polyclonal Antibody	A58370	Human	ELISA, WB
C4A Polyclonal Antibody	A69415	Human	ELISA, WB, IHC
C4B Polyclonal Antibody	A56158	Human	ELISA, IHC
C4BPA Polyclonal Antibody	A62178	Human	ELISA, IHC
C4orf3 Polyclonal Antibody	A56819	Human	ELISA, IHC, IF
C4orf3 Polyclonal Antibody	A66426	Human	ELISA, WB
C4orf33 Polyclonal Antibody	A66338	Human, Mouse	ELISA, WB
C4orf46 Polyclonal Antibody	A65074	Human	ELISA, IHC
C5 Polyclonal Antibody	A53695	Pig, Human	ELISA, IHC
C5 Polyclonal Antibody	A58374	Human	ELISA, IHC
C5orf22 Polyclonal Antibody	A58378	Human, Mouse	ELISA, WB, IHC
C5orf24 Polyclonal Antibody	A65898	Human, Mouse	ELISA, WB, IHC
C5orf28 Polyclonal Antibody	A64618	Human	ELISA, IF
C5orf45 Polyclonal Antibody	A65470	Human, Mouse	ELISA, WB
C5orf51 Polyclonal Antibody	A62182	Human	ELISA, IHC, IF
C6orf136 Polyclonal Antibody	A65162	Human	ELISA
C6orf141 Polyclonal Antibody	A65258	Human	ELISA, IHC
C6orf203 Polyclonal Antibody	A58382	Human	ELISA, IHC
C6orf89 Polyclonal Antibody	A65714	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
C7orf33 Polyclonal Antibody	A58386	Human	ELISA
C7orf34 Polyclonal Antibody	A66602	Human	ELISA, WB, IHC
C7orf50 Polyclonal Antibody	A51273	Human	ELISA, WB
C7orf61 Polyclonal Antibody	A58390	Human	ELISA, IHC
C7orf62 Polyclonal Antibody	A58394	Human	ELISA, WB, IHC
C8G Polyclonal Antibody	A57261	Human	ELISA, WB
C8orf22 Polyclonal Antibody	A66666	Human	ELISA, IHC
C8orf31 Polyclonal Antibody	A66758	Human	ELISA
C8orf34 Polyclonal Antibody	A65002	Human, Mouse	ELISA, WB, IHC
C8orf48 Polyclonal Antibody	A58398	Human, Mouse	ELISA, WB, IHC
C8orf59 Polyclonal Antibody	A66478	Human	ELISA
C8orf74 Polyclonal Antibody	A65686	Human	ELISA, IHC
C9orf117 Polyclonal Antibody	A65062	Human	ELISA, IHC
C9orf16 Polyclonal Antibody	A58402	Human	ELISA, IHC
C9orf163 Polyclonal Antibody	A56984	Human	ELISA, IHC
C9orf43 Polyclonal Antibody	A66230	Human	ELISA, IHC
C9orf57 Polyclonal Antibody	A65298	Human	ELISA, IHC
C9orf64 Polyclonal Antibody	A65166	Human	ELISA, WB, IHC
C9orf72 Polyclonal Antibody	A66710	Human	ELISA, IHC, IF
Ca1 Polyclonal Antibody	A53734	Rat	ELISA, WB
Ca1 Polyclonal Antibody	A54816	Mouse, Rat	ELISA, WB
CA1 Polyclonal Antibody	A55202	Human, Mouse, Rat	ELISA, WB, IHC, IF
CA11 Polyclonal Antibody	A58410	Human	ELISA, IHC
CA12 Polyclonal Antibody	A55671	Human, Mouse	ELISA, WB, IHC
CA12 Polyclonal Antibody	A55882	Human, Mouse	ELISA, WB
CA2 Polyclonal Antibody	A55113	Human, Mouse, Rat	ELISA, WB, IHC, IF
CA3 Polyclonal Antibody	A62194	Human, Mouse	ELISA, WB, IHC
CA5B Polyclonal Antibody	A51417	Human, Mouse	ELISA, WB, IHC
CAB39L Polyclonal Antibody	A58414	Human	ELISA, WB
CABP4 Polyclonal Antibody	A69419	Human	ELISA, IHC, IF
CABP5 Polyclonal Antibody	A67294	Human, Mouse	ELISA, WB
CACNA2D1 Polyclonal Antibody	A56306	Human	ELISA, IHC
CACNA2D4 Polyclonal Antibody	A65974	Human	ELISA, IHC
CACNB2 Polyclonal Antibody	A69031	Human	ELISA, IHC, IF
cagA Polyclonal Antibody	A52066	Helicobacter pylori	ELISA
cagA Polyclonal Antibody	A57852	Helicobacter pylori	ELISA
cagA Polyclonal Antibody	A64204	Helicobacter pylori	ELISA
CALCA Polyclonal Antibody	A52257	Human	ELISA
CALHM3 Polyclonal Antibody	A65770	Human, Mouse	ELISA, WB, IHC, IF
CALM1 Polyclonal Antibody	A51453	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
CALM1 Polyclonal Antibody	A55511	Human, Mouse	ELISA, WB
CALM1 Polyclonal Antibody	A55759	Human	ELISA
CALM1 Polyclonal Antibody	A58418	Human	ELISA, IF
CALML4 Polyclonal Antibody	A51077	Human	ELISA, IHC
CALML5 Polyclonal Antibody	A51313	Human	ELISA, WB, IHC
CALN1 Polyclonal Antibody	A51297	Human, Mouse	ELISA, WB
CALR Polyclonal Antibody	A50640	Human	ELISA, WB, IHC
CALR Polyclonal Antibody	A52608	Human	ELISA, WB, IHC
CALR Polyclonal Antibody	A69423	Pig	ELISA
Camk2n1 Polyclonal Antibody	A58422	Mouse	ELISA
Canine Serum Albumin Polyclonal Antibody	A57564	Canine	ELISA
CANT1 Polyclonal Antibody	A55890	Human	ELISA
CAP2 Polyclonal Antibody	A62198	Human, Mouse	ELISA, WB, IHC
CAPN1 Polyclonal Antibody	A57493	Human	ELISA, WB, IHC
CAPN2 Polyclonal Antibody	A52892	Human, Mouse	ELISA, WB
CAPN2 Polyclonal Antibody	A57049	Pig	ELISA
CAPN6 Polyclonal Antibody	A68238	Human	ELISA, WB, IHC
CAPNS1 Polyclonal Antibody	A56675	Human	ELISA
Caprin1 Polyclonal Antibody	A56923	Rat	ELISA
CAPS2 Polyclonal Antibody	A69171	Human	ELISA, WB, IHC
Capsid Protein VP2 Polyclonal Antibody	A57197	Human parvovirus B19	ELISA
CAPZA2 Polyclonal Antibody	A50093	Human	ELISA, IHC
CARD19 Polyclonal Antibody	A58406	Human	ELISA, IHC
CARHSP1 Polyclonal Antibody	A58426	Human	ELISA
CARS Polyclonal Antibody	A62202	Human, Mouse	ELISA, WB, IHC
Casein Polyclonal Antibody	A57806	Human	ELISA
CASK Polyclonal Antibody	A58430	Human	ELISA, IHC
CASP1 Polyclonal Antibody	A51865	Human	ELISA, IHC
CASP1 Polyclonal Antibody	A64084	Human	ELISA, IHC
CASP10 Polyclonal Antibody	A66574	Human	ELISA
Casp12 Polyclonal Antibody	A50097	Mouse, Human	ELISA, WB
CASP2 Polyclonal Antibody	A57285	Human	ELISA, WB
CASP3 Polyclonal Antibody	A52905	Human, Mouse	ELISA, WB, IHC
CASP7 Polyclonal Antibody	A54193	Human	ELISA, WB
CASP8 Polyclonal Antibody	A53076	Human	ELISA, WB, IHC, IF
CASQ1 Polyclonal Antibody	A62206	Human, Mouse	ELISA, WB, IHC
CASR Polyclonal Antibody	A53024	Human	ELISA, IF
CAST Polyclonal Antibody	A54656	Human	ELISA, WB, IHC
Cathepsin D Polyclonal Antibody	A50157	Cricetulus griseus	ELISA
CATSPER2 Polyclonal Antibody	A58434	Human, Mouse	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
CATSPER4 Polyclonal Antibody	A65454	Human	ELISA, IHC, IF
CBFB Polyclonal Antibody	A58438	Human	ELISA
Cbs Polyclonal Antibody	A66894	Mouse, Human	ELISA, IF
CBWD2 Polyclonal Antibody	A66118	Human, Mouse	ELISA, WB, IHC, IF
CBX2 Polyclonal Antibody	A58442	Human, Mouse	ELISA, WB, IHC
CBX5 Polyclonal Antibody	A50101	Human	ELISA, WB, IHC
CCDC103 Polyclonal Antibody	A66098	Human, Mouse	ELISA, WB
CCDC106 Polyclonal Antibody	A58450	Human, Mouse	ELISA, WB, IHC
CCDC108 Polyclonal Antibody	A58454	Human	ELISA, WB, IHC, IF
CCDC109B Polyclonal Antibody	A58458	Human	ELISA, IHC, IF
CCDC112 Polyclonal Antibody	A58462	Human	ELISA, WB, IHC, IF
CCDC113 Polyclonal Antibody	A67682	Human	ELISA, IHC
CCDC116 Polyclonal Antibody	A58466	Human	ELISA, WB, IHC
CCDC127 Polyclonal Antibody	A58470	Human, Mouse	ELISA, WB
CCDC134 Polyclonal Antibody	A51245	Human	ELISA, IHC
CCDC14 Polyclonal Antibody	A50953	Human	ELISA, WB, IHC
CCDC150 Polyclonal Antibody	A66506	Human	ELISA, IHC
CCDC155 Polyclonal Antibody	A66194	Human	ELISA
CCDC173 Polyclonal Antibody	A64630	Human	ELISA
CCDC181 Polyclonal Antibody	A58474	Human	ELISA, IHC
CCDC184 Polyclonal Antibody	A65038	Human	ELISA, IHC
CCDC24 Polyclonal Antibody	A58478	Human	ELISA, WB, IHC
CCDC36 Polyclonal Antibody	A51037	Human	ELISA, IHC
CCDC42 Polyclonal Antibody	A58482	Human, Mouse	ELISA, WB, IHC
CCDC53 Polyclonal Antibody	A58486	Human	ELISA
CCDC57 Polyclonal Antibody	A58490	Human	ELISA, WB, IHC
CCDC59 Polyclonal Antibody	A67766	Human	ELISA, IHC
CCDC66 Polyclonal Antibody	A62210	Human	ELISA, IHC
CCDC68 Polyclonal Antibody	A58494	Human	ELISA, IHC
CCDC79 Polyclonal Antibody	A66382	Human	ELISA, IHC
CCDC82 Polyclonal Antibody	A58498	Human	ELISA
CCDC84 Polyclonal Antibody	A65846	Human	ELISA, IHC
CCDC87 Polyclonal Antibody	A67614	Human	ELISA
CCDC89 Polyclonal Antibody	A66210	Human, Mouse	ELISA, WB, IHC
CCDC91 Polyclonal Antibody	A58502	Human	ELISA, IHC
CCDC94 Polyclonal Antibody	A67346	Human	ELISA, IHC
CCDC97 Polyclonal Antibody	A58506	Human	ELISA, IHC
Ccl12 Polyclonal Antibody	A53181	Mouse	ELISA
CCL14 Polyclonal Antibody	A55419	Human	ELISA, IHC
CCL16 Polyclonal Antibody	A55423	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
CCL2 Polyclonal Antibody	A51548	Human	ELISA, WB, IHC, IF
Ccl2 Polyclonal Antibody	A51954	Mouse, Human	ELISA, IF
CCL2 Polyclonal Antibody	A52218	Human	ELISA, IHC, IF
CCL21 Polyclonal Antibody	A53746	Human	ELISA
CCL22 Polyclonal Antibody	A51445	Human	ELISA, WB
CCL24 Polyclonal Antibody	A54017	Human	ELISA, IHC
Ccl25 Polyclonal Antibody	A53085	Mouse	ELISA
CCL4 Polyclonal Antibody	A53170	Mouse, Human	ELISA, IHC
CCL5 Polyclonal Antibody	A51637	Human	ELISA, WB
CCL5 Polyclonal Antibody	A52420	Human	ELISA, WB
CCL7 Polyclonal Antibody	A51693	Human	ELISA, WB
CCL8 Polyclonal Antibody	A54970	Human	ELISA, IHC
Ccl9 Polyclonal Antibody	A53404	Mouse	ELISA
CCND1 Polyclonal Antibody	A54316	Human	ELISA
CCNJL Polyclonal Antibody	A58510	Human, Mouse	ELISA, WB, IHC
CCNY Polyclonal Antibody	A70111	Human	ELISA, WB, IHC, IF
CCR1 Polyclonal Antibody	A62214	Human	ELISA, IHC, IF
CCR4 Polyclonal Antibody	A62218	Human	ELISA, IF
Ccr5 Polyclonal Antibody	A58514	Mouse	ELISA
CCR6 Polyclonal Antibody	A50105	Human	ELISA, WB
CCS Polyclonal Antibody	A68475	Human	ELISA, IHC
CCT6B Polyclonal Antibody	A58518	Human	ELISA, IHC
CD14 Polyclonal Antibody	A64104	Human	ELISA, IHC
CD160 Polyclonal Antibody	A50109	Human	ELISA, IHC
Cd163 Polyclonal Antibody	A53832	Mouse, Human	ELISA, WB
CD177 Polyclonal Antibody	A66950	Human, Mouse	ELISA, WB, IHC, IF
CD1E Polyclonal Antibody	A68479	Human	ELISA, IHC
CD2 Polyclonal Antibody	A57021	Human	ELISA, IHC
CD27 Polyclonal Antibody	A50648	Human	ELISA
CD300LB Polyclonal Antibody	A62222	Human	ELISA
CD300LF Polyclonal Antibody	A57477	Human	ELISA, IHC
CD300LG Polyclonal Antibody	A68975	Human, Rat	ELISA, WB, IF
CD302 Polyclonal Antibody	A66166	Human	ELISA
CD3E Polyclonal Antibody	A51946	Human	ELISA, WB
CD40LG Polyclonal Antibody	A50624	Human	ELISA
CD46 Polyclonal Antibody	A50113	Human	ELISA, WB, IHC
Cd46 Polyclonal Antibody	A53651	Mouse, Human	ELISA, WB, IHC
CD59 Polyclonal Antibody	A53261	Human	ELISA, IHC
CD59 Polyclonal Antibody	A53551	Human, Mouse	ELISA, WB, IHC
Cd63 Polyclonal Antibody	A53109	Mouse, Human	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
Cd68 Polyclonal Antibody	A56178	Mouse	ELISA
CD7 Polyclonal Antibody	A55667	Human	ELISA, IHC
Cd74 Polyclonal Antibody	A54780	Mouse	ELISA, WB
Cd74 Polyclonal Antibody	A55559	Rat, Mouse	ELISA, WB
Cd81 Polyclonal Antibody	A57217	Mouse	ELISA
Cd82 Polyclonal Antibody	A55487	Mouse, Human	ELISA, WB
CD86 Polyclonal Antibody	A50117	Human	ELISA
CD8A Polyclonal Antibody	A62226	Human	ELISA, IHC
CD99L2 Polyclonal Antibody	A69083	Human	ELISA, IHC, IF
CDA Polyclonal Antibody	A52901	Human	ELISA, WB, IHC, IF
CDADC1 Polyclonal Antibody	A67542	Human	ELISA, IHC
CDC25C Polyclonal Antibody	A58522	Human	ELISA, IHC
CDC26 Polyclonal Antibody	A66406	Human	ELISA, IHC
CDC34 Polyclonal Antibody	A62230	Human	ELISA, IHC
CDC42 Polyclonal Antibody	A62234	Human	ELISA, WB
CDC42BPB Polyclonal Antibody	A68787	Human	ELISA, IHC, IF
CDC42EP2 Polyclonal Antibody	A62238	Human	ELISA, IHC
CDCA7L Polyclonal Antibody	A70139	Human	ELISA, IHC
CDH12 Polyclonal Antibody	A52170	Human	ELISA, WB, IHC, IF
CDH26 Polyclonal Antibody	A66170	Human	ELISA
CDH3 Polyclonal Antibody	A50660	Human	ELISA, IHC, IF
CDH7 Polyclonal Antibody	A69035	Human	ELISA, WB, IHC, IF
CDH9 Polyclonal Antibody	A69319	Human, Mouse	ELISA, WB, IHC, IF
CDK11A Polyclonal Antibody	A54009	Human	ELISA, WB, IHC, IF
CDK12 Polyclonal Antibody	A70299	Human	ELISA, WB
CDK2 Polyclonal Antibody	A62242	Human	ELISA, WB, IHC, IF, IP
CDK2AP1 Polyclonal Antibody	A53256	Human	ELISA, IHC
CDK2AP2 Polyclonal Antibody	A62246	Human	ELISA, IHC
CDK4 Polyclonal Antibody	A52392	Human	ELISA, IHC
CDK5 Polyclonal Antibody	A50121	Human, Mouse	ELISA, WB, IHC, IF
CDKB1-2 Polyclonal Antibody	A64874	Arabidopsis thaliana	ELISA
CDKN1A Polyclonal Antibody	A53012	Human	ELISA
CDKN2AIPNL Polyclonal Antibody	A53293	Human, Mouse	ELISA, IHC
CDKN2B Polyclonal Antibody	A62250	Human	ELISA, IHC, IF
CDPF1 Polyclonal Antibody	A65682	Human	ELISA
CDR2 Polyclonal Antibody	A52560	Human, Mouse	ELISA, WB, IHC
CDR2L Polyclonal Antibody	A69427	Human	ELISA, WB, IHC
CEACAM1 Polyclonal Antibody	A58526	Human	ELISA, WB
CEACAM16 Polyclonal Antibody	A69431	Human	ELISA, IHC
CEACAM7 Polyclonal Antibody	A62254	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
CEL Polyclonal Antibody	A69823	Human	ELISA, IHC
Cela2a Polyclonal Antibody	A54384	Mouse	ELISA, WB
CELA3A Polyclonal Antibody	A56968	Human, Mouse	ELISA, WB
CENPB Polyclonal Antibody	A50053	Human	ELISA, IHC
CENPJ Polyclonal Antibody	A69435	Human	ELISA, WB, IF
CEP126 Polyclonal Antibody	A55794	Human	ELISA, IHC
CEP135 Polyclonal Antibody	A68927	Human	ELISA, IF
CEP250 Polyclonal Antibody	A70235	Human	ELISA, IF
CEP63 Polyclonal Antibody	A51101	Human, Mouse	ELISA, WB, IHC
CEP68 Polyclonal Antibody	A65738	Human	ELISA
CEP76 Polyclonal Antibody	A58530	Human	ELISA, WB, IHC
CEP83 Polyclonal Antibody	A56787	Human	ELISA
CERS3 Polyclonal Antibody	A70003	Human, Mouse	ELISA, WB
CES1 Polyclonal Antibody	A62258	Human, Mouse	ELISA, WB, IHC, IF
Ces1c Polyclonal Antibody	A53546	Human, Mouse, Rat	ELISA, WB, IHC
CETN2 Polyclonal Antibody	A54898	Human	ELISA
cfaE Polyclonal Antibody	A64322	Escherichia coli	ELISA
CFAP100 Polyclonal Antibody	A64978	Human	ELISA, IHC, IF
CFAP52 Polyclonal Antibody	A66614	Human	ELISA, WB, IHC
CFAP61 Polyclonal Antibody	A58534	Human	ELISA, IHC
CFB Polyclonal Antibody	A55575	Human	ELISA, WB
CFHR2 Polyclonal Antibody	A58538	Human	ELISA, WB, IHC
CFLAR Polyclonal Antibody	A58542	Human	ELISA, WB, IHC
CGGBP1 Polyclonal Antibody	A68018	Human	ELISA, IHC
CHAC1 Polyclonal Antibody	A58546	Human, Mouse	ELISA, WB, IHC
CHAD Polyclonal Antibody	A58550	Human	ELISA, IHC
CHCHD4 Polyclonal Antibody	A56314	Human	ELISA, IHC
CHEK2 Polyclonal Antibody	A62262	Human	ELISA, WB
CHGA Polyclonal Antibody	A54930	Human	ELISA, IHC
CHI3L2 Polyclonal Antibody	A55814	Human	ELISA, IHC
Chicken Serum Albumin Polyclonal Antibody	A57529	Chicken	ELISA
CHID1 Polyclonal Antibody	A58554	Human	ELISA, WB, IHC
CHM Polyclonal Antibody	A62266	Human	ELISA, IHC
CHMP1A Polyclonal Antibody	A69259	Human	ELISA, IF
CHMP2A Polyclonal Antibody	A51805	Human	ELISA, WB, IHC
CHMP5 Polyclonal Antibody	A57253	Human	ELISA, WB
CHMP6 Polyclonal Antibody	A58558	Human, Rat	ELISA, WB, IHC
CHMP7 Polyclonal Antibody	A58562	Human	ELISA, IHC
CHN2 Polyclonal Antibody	A68483	Human	ELISA, IHC
choD Polyclonal Antibody	A55778	Mycobacterium tuberculosis	ELISA

Product Name	Cat. #	Species Reactivity	Applications
choD Polyclonal Antibody	A64196	Mycobacterium tuberculosis	ELISA
Chrm1 Polyclonal Antibody	A53520	Rat	ELISA
CHRM3 Polyclonal Antibody	A57145	Human	ELISA
CHRNA1 Polyclonal Antibody	A53586	Human, Mouse, Rat	ELISA, WB, IHC
CHRNA1 Polyclonal Antibody	A54824	Torpedo californica	ELISA, WB
Chrna1 Polyclonal Antibody	A55463	Mouse, Human	ELISA, WB
CHRN2 Polyclonal Antibody	A68487	Human, Rat	ELISA, WB, IF
CHRN3 Polyclonal Antibody	A68491	Human, Mouse	ELISA, WB, IHC, IF
CHS1 Polyclonal Antibody	A54633	Baker's yeast	ELISA
CHST10 Polyclonal Antibody	A58566	Human	ELISA
CHST11 Polyclonal Antibody	A70291	Human	ELISA, IHC
CIAPIN1 Polyclonal Antibody	A57093	Human	ELISA
CIART Polyclonal Antibody	A58570	Human	ELISA, IHC
CINP Polyclonal Antibody	A51229	Human	ELISA
CIPC Polyclonal Antibody	A67674	Human, Mouse	ELISA, WB, IHC
CISD2 Polyclonal Antibody	A66626	Human, Mouse	ELISA, WB, IHC, IF
CKAP5 Polyclonal Antibody	A69859	Human, Rat	ELISA, WB, IHC, IF
CKB Polyclonal Antibody	A50382	Human, Mouse	ELISA, WB, IHC, IF
CKM Polyclonal Antibody	A52021	Human, Mouse	ELISA, WB, IHC
CKMT1A Polyclonal Antibody	A52297	Human	ELISA, IHC, IF
CKS2 Polyclonal Antibody	A62270	Human	ELISA, IHC
CLASP2 Polyclonal Antibody	A58574	Human	ELISA, IHC
CLC Polyclonal Antibody	A69795	Human	ELISA, IHC, IF
CLCN2 Polyclonal Antibody	A62274	Human	ELISA, IF
CLCNKB Polyclonal Antibody	A69439	Human, Rat	ELISA, WB, IHC, IF
CLDN12 Polyclonal Antibody	A62278	Human	ELISA, IHC
CLDN15 Polyclonal Antibody	A50125	Human	ELISA, IHC, IF
CLDN20 Polyclonal Antibody	A62282	Human	ELISA, WB, IF
CLEC18A Polyclonal Antibody	A54475	Human, Mouse	ELISA, WB, IHC
CLEC1A Polyclonal Antibody	A58578	Human, Mouse	ELISA, WB, IHC
CLEC2D Polyclonal Antibody	A67778	Human	ELISA
CLEC3A Polyclonal Antibody	A62286	Human, Mouse	ELISA, WB, IHC
CLEC4C Polyclonal Antibody	A58582	Human	ELISA, IHC
CLEC4D Polyclonal Antibody	A57325	Human	ELISA, IHC
CLEC4E Polyclonal Antibody	A58586	Human, Mouse	ELISA, WB
clfB Polyclonal Antibody	A65546	Staphylococcus aureus	ELISA
CLGN Polyclonal Antibody	A62290	Human, Mouse	ELISA, WB, IHC
CLIC2 Polyclonal Antibody	A62294	Human	ELISA, IHC, IF
CLN5 Polyclonal Antibody	A68495	Human	ELISA, IF
CLPB Polyclonal Antibody	A58590	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
CLPP Polyclonal Antibody	A68851	Human	ELISA, IHC
CLPS Polyclonal Antibody	A69443	Human	ELISA, IHC
clpS Polyclonal Antibody	A52134	Escherichia coli	ELISA, WB
CLPTM1L Polyclonal Antibody	A70143	Human	ELISA, WB, IHC, IF
CLRN2 Polyclonal Antibody	A62298	Human	ELISA, IHC
CLRN3 Polyclonal Antibody	A66218	Human	ELISA, IF
CLSTN2 Polyclonal Antibody	A67266	Human	ELISA, IHC, IF
CLTB Polyclonal Antibody	A55735	Human	ELISA, IHC
CLTCL1 Polyclonal Antibody	A55178	Human	ELISA, WB, IHC
CLU Polyclonal Antibody	A52202	Human	ELISA, WB, IHC
CLUAP1 Polyclonal Antibody	A58594	Human	ELISA, IHC
CLVS2 Polyclonal Antibody	A56471	Human, Mouse	ELISA, WB
CLYBL Polyclonal Antibody	A66938	Human, Mouse	ELISA, WB, IHC
Cma1 Polyclonal Antibody	A55523	Mouse	ELISA
CMA1 Polyclonal Antibody	A62302	Human, Mouse	ELISA, WB
CMBL Polyclonal Antibody	A58598	Human, Mouse	ELISA, WB, IHC, IF
CMC1 Polyclonal Antibody	A62306	Human	ELISA, IHC
CMC4 Polyclonal Antibody	A64410	Human	ELISA, IHC
cmk Polyclonal Antibody	A50845	Escherichia coli	ELISA
CMTM2 Polyclonal Antibody	A58602	Human	ELISA, IF
CMTM5 Polyclonal Antibody	A66694	Human	ELISA, IHC
CMTM8 Polyclonal Antibody	A66054	Human	ELISA, IF
CMTR2 Polyclonal Antibody	A58606	Human	ELISA, IHC, IF
CNBP Polyclonal Antibody	A50129	Human	ELISA, IHC
CNDP2 Polyclonal Antibody	A54304	Human	ELISA, IHC
CNGA4 Polyclonal Antibody	A66018	Human	ELISA, IHC
CNIH2 Polyclonal Antibody	A65590	Human, Mouse	ELISA, WB, IHC, IF
CNIH3 Polyclonal Antibody	A66870	Human	ELISA, IHC, IF
CNIH4 Polyclonal Antibody	A58610	Human	ELISA, WB, IHC, IF
CNKS2 Polyclonal Antibody	A57125	Human	ELISA, IHC
CNNM1 Polyclonal Antibody	A67602	Human	ELISA, IF
CNOT2 Polyclonal Antibody	A68262	Human	ELISA, WB
CNPY2 Polyclonal Antibody	A53985	Human	ELISA, WB, IHC
CNPY4 Polyclonal Antibody	A58618	Human	ELISA, IHC
CNTF Polyclonal Antibody	A52282	Human	ELISA, IHC
CNTFR Polyclonal Antibody	A50133	Human	ELISA, WB, IHC
CNTNAP1 Polyclonal Antibody	A50137	Human	ELISA, IHC
COA1 Polyclonal Antibody	A62186	Human	ELISA, IHC, IF
COA6 Polyclonal Antibody	A65050	Human	ELISA, IHC
COBLL1 Polyclonal Antibody	A58622	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
COG3 Polyclonal Antibody	A66598	Human	ELISA, IHC
COL12A1 Polyclonal Antibody	A55162	Chicken	ELISA
COL12A1 Polyclonal Antibody	A67142	Human	ELISA, IHC
COL17A1 Polyclonal Antibody	A52130	Human	ELISA, IHC
Col1a1 Polyclonal Antibody	A53639	Rat, Human	ELISA, IHC
COL20A1 Polyclonal Antibody	A58626	Human	ELISA, IHC
COL2A1 Polyclonal Antibody	A54826	Human	ELISA, IHC
COL3A1 Polyclonal Antibody	A54145	Human	ELISA
COL4A1 Polyclonal Antibody	A54149	Human	ELISA, IHC, IF
COL4A1 Polyclonal Antibody	A55058	Human	ELISA, IHC
COL4A2 Polyclonal Antibody	A54161	Human	ELISA, IHC
COL4A3 Polyclonal Antibody	A57882	Human	ELISA, IHC
COL4A3BP Polyclonal Antibody	A68194	Human, Mouse	ELISA, WB, IHC, IF
COL6A6 Polyclonal Antibody	A50865	Human	ELISA, IHC
COL6A6 Polyclonal Antibody	A57485	Human	ELISA, IHC
COL8A1 Polyclonal Antibody	A62310	Human, Mouse	ELISA, WB, IF
COLEC10 Polyclonal Antibody	A56980	Human, Mouse	ELISA, WB, IHC
COLGALT2 Polyclonal Antibody	A58630	Human, Mouse	ELISA, WB, IHC
COMMD2 Polyclonal Antibody	A65926	Human	ELISA, IHC
COMMD4 Polyclonal Antibody	A51921	Human	ELISA, IHC
COMMD6 Polyclonal Antibody	A65854	Human	ELISA, IHC
COMMD9 Polyclonal Antibody	A58634	Human	ELISA, IHC
Conglutin Polyclonal Antibody	A65218	Peanut	ELISA
COPB1 Polyclonal Antibody	A58638	Human, Mouse	ELISA, WB, IHC
COPE Polyclonal Antibody	A64056	Human	ELISA, WB, IHC
COPG2 Polyclonal Antibody	A70215	Human	ELISA, IHC, IF
COPS7A Polyclonal Antibody	A56815	Human, Mouse	ELISA, WB, IHC
COPS7B Polyclonal Antibody	A67514	Human, Mouse	ELISA, WB, IHC, IF
COPS8 Polyclonal Antibody	A69163	Human	ELISA, IHC
COPZ1 Polyclonal Antibody	A62314	Human, Mouse	ELISA, WB, IHC
COQ4 Polyclonal Antibody	A68182	Human	ELISA, IHC
CORO7 Polyclonal Antibody	A62318	Human	ELISA
COTL1 Polyclonal Antibody	A55651	Human, Mouse	ELISA, WB, IHC
COX11 Polyclonal Antibody	A58642	Human	ELISA, IHC, IF
COX17 Polyclonal Antibody	A64770	Human	ELISA, WB, IHC
COX19 Polyclonal Antibody	A65026	Human	ELISA, IF
COX4I1 Polyclonal Antibody	A55571	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC
COX5A Polyclonal Antibody	A52616	Human	ELISA, WB, IHC
COX5B Polyclonal Antibody	A51597	Human	ELISA, WB, IF
COX6A1 Polyclonal Antibody	A52685	Human	ELISA, WB, IHC, IF



Product Name	Cat. #	Species Reactivity	Applications
COX6B1 Polyclonal Antibody	A51609	Human	ELISA, IF
CP Polyclonal Antibody	A57640	Human	ELISA, IHC
CPA2 Polyclonal Antibody	A58646	Human	ELISA, WB, IHC
CPA6 Polyclonal Antibody	A50141	Human	ELISA
CPE Polyclonal Antibody	A69447	Human	ELISA, WB
CPEB1 Polyclonal Antibody	A51233	Human	ELISA
CPEB4 Polyclonal Antibody	A58650	Human	ELISA, WB
CPED1 Polyclonal Antibody	A58654	Human	ELISA, IHC
cpg2 Polyclonal Antibody	A54644	Pseudomonas sp	ELISA, WB
CPLX1 Polyclonal Antibody	A62322	Human	ELISA, IHC, IF
CPLX3 Polyclonal Antibody	A58658	Human	ELISA
CPLX4 Polyclonal Antibody	A65750	Human	ELISA
CPNE2 Polyclonal Antibody	A66590	Human	ELISA, IHC
CPNE4 Polyclonal Antibody	A66906	Human, Mouse, Rat	ELISA, WB, IHC, IF
CPNE9 Polyclonal Antibody	A58662	Human	ELISA, IHC
CPOX Polyclonal Antibody	A68499	Human	ELISA, WB, IF
CPS1 Polyclonal Antibody	A69451	Human	ELISA, WB, IHC, IF
CPVL Polyclonal Antibody	A55934	Human	ELISA, IHC
CPXCR1 Polyclonal Antibody	A66334	Human, Mouse	ELISA, WB, IHC
Cr2 Polyclonal Antibody	A62326	Mouse	ELISA
CRABP1 Polyclonal Antibody	A55078	Human, Mouse	ELISA, WB, IHC
CRABP2 Polyclonal Antibody	A62330	Human	ELISA, WB, IHC
CRADD Polyclonal Antibody	A55663	Human	ELISA, WB
CREB1 Polyclonal Antibody	A52845	Human, Mouse	ELISA, WB, IHC
creN7 Polyclonal Antibody	A55906	Sulfolobus islandicus	ELISA
CRHR1 Polyclonal Antibody	A53655	Human, Mouse	ELISA, WB, IHC, IF
CRHR1 Polyclonal Antibody	A55411	Human	ELISA
CRIP2 Polyclonal Antibody	A57369	Human	ELISA
CRIPT Polyclonal Antibody	A62334	Human	ELISA
CRISPLD1 Polyclonal Antibody	A67826	Human	ELISA, IHC
CRLF1 Polyclonal Antibody	A69455	Human	ELISA, IHC
CRMP1 Polyclonal Antibody	A56927	Human, Mouse, Rat	ELISA, WB
CRNN Polyclonal Antibody	A58666	Human, Mouse, Rat	ELISA, WB
crp Polyclonal Antibody	A55435	Escherichia coli	ELISA
CRP Polyclonal Antibody	A52500	Human	ELISA, WB, IHC
cry1Ab Polyclonal Antibody	A55098	Bacillus thuringiensis kurstaki	ELISA
cry1Ab Polyclonal Antibody	A55527	Bacillus thuringiensis kurstaki	ELISA
cry1Ac Polyclonal Antibody	A54664	Bacillus thuringiensis kurstaki	ELISA, WB
cry1Ac Polyclonal Antibody	A55038	Bacillus thuringiensis	ELISA
cry1Fb Polyclonal Antibody	A54552	Bacillus thuringiensis morrisoni	ELISA

Product Name	Cat. #	Species Reactivity	Applications
cry3Aa Polyclonal Antibody	A54551	Bacillus thuringiensis tenebrionis	ELISA
CRYAA Polyclonal Antibody	A68503	Human	ELISA, IF
CRYAB Polyclonal Antibody	A55834	Human, Mouse, Rat	ELISA, WB, IHC
CRYBA1 Polyclonal Antibody	A62338	Human, Mouse	ELISA, WB
CRYBA4 Polyclonal Antibody	A62342	Human, Mouse	ELISA, WB, IHC
CRYGN Polyclonal Antibody	A66670	Human	ELISA, IHC
CRYM Polyclonal Antibody	A68891	Human	ELISA, IHC, IF
CRYZL1 Polyclonal Antibody	A58670	Human	ELISA, WB
CS Polyclonal Antibody	A52776	Human, Mouse	ELISA, WB, IHC
Csad Polyclonal Antibody	A58674	Mouse	ELISA, WB
CSAG1 Polyclonal Antibody	A62346	Human	ELISA, IHC
Csf1 Polyclonal Antibody	A51905	Mouse	ELISA
Csf1 Polyclonal Antibody	A52920	Mouse, Human	ELISA, IHC
CSF1 Polyclonal Antibody	A56667	Human	ELISA, IHC, IF
CSF1R Polyclonal Antibody	A54832	Felis silvestris catus	ELISA
CSF2 Polyclonal Antibody	A55375	Human	ELISA, IHC
CSF2RA Polyclonal Antibody	A56976	Human	ELISA, IHC
CSF3 Polyclonal Antibody	A52476	Human	ELISA, IHC
CSGALNACT2 Polyclonal Antibody	A58678	Human	ELISA, WB, IHC
CSH1 Polyclonal Antibody	A57619	Human	ELISA, WB
CSHL1 Polyclonal Antibody	A50145	Human, Mouse	ELISA, WB, IHC
CSMD2 Polyclonal Antibody	A65826	Human	ELISA, IHC
CSMD3 Polyclonal Antibody	A69991	Human	ELISA, IHC, IF
CSN2 Polyclonal Antibody	A69459	Human	ELISA, IHC
CSNK1A1 Polyclonal Antibody	A54428	Human	ELISA, IHC
CSNK1D Polyclonal Antibody	A58682	Human	ELISA, IHC
CSNK1E Polyclonal Antibody	A52740	Human	ELISA, WB
CSNK1E Polyclonal Antibody	A56246	Human	ELISA, IHC, IF
CSNK2A1 Polyclonal Antibody	A57361	Human	ELISA, WB, IHC, IF
CSNK2A2 Polyclonal Antibody	A54748	Human	ELISA, WB, IHC
CSNK2B Polyclonal Antibody	A51553	Human	ELISA, IHC
CSRP1 Polyclonal Antibody	A51966	Human	ELISA, IHC
cssB Polyclonal Antibody	A50804	Escherichia coli	ELISA
CST11 Polyclonal Antibody	A67938	Human	ELISA, IHC
CST3 Polyclonal Antibody	A53373	Human	ELISA, WB, IHC
CST9L Polyclonal Antibody	A67186	Human	ELISA, IHC
CSTA Polyclonal Antibody	A50149	Human	ELISA, WB, IHC
CSTB Polyclonal Antibody	A52620	Human	ELISA, WB, IHC
CTDNEP1 Polyclonal Antibody	A54013	Human	ELISA, IHC
CTGF Polyclonal Antibody	A52384	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
CTHRC1 Polyclonal Antibody	A51065	Human	ELISA, IHC
CTNNA1 Polyclonal Antibody	A58686	Human, Mouse	ELISA, WB
CTNNB1 Polyclonal Antibody	A51457	Human, Mouse, Zebrafish	ELISA, WB, IHC
CTPS1 Polyclonal Antibody	A50153	Human	ELISA, WB, IHC
CTPS2 Polyclonal Antibody	A67302	Human	ELISA, WB, IHC
CTRB1 Polyclonal Antibody	A58690	Human, Mouse	ELISA, WB, IHC
CTSA Polyclonal Antibody	A62350	Human	ELISA, IHC, IF
CTSB Polyclonal Antibody	A51869	Human, Mouse	ELISA, WB, IHC, IF
CTSD Polyclonal Antibody	A52324	Human	ELISA
CTSF Polyclonal Antibody	A52114	Human	ELISA
Ctsg Polyclonal Antibody	A57421	Mouse	ELISA
CTSK Polyclonal Antibody	A52840	Human	ELISA
Ctsk Polyclonal Antibody	A53433	Rat, Human	ELISA, IHC
CTSL Polyclonal Antibody	A50161	Chlorocebus aethiops	ELISA, WB
CTSS Polyclonal Antibody	A53500	Human	ELISA, IHC
Ctss Polyclonal Antibody	A56507	Mouse	ELISA
CUEDC2 Polyclonal Antibody	A56763	Human	ELISA
CUTC Polyclonal Antibody	A70243	Human	ELISA, IHC, IF
CUX1 Polyclonal Antibody	A69855	Human, Mouse, Rat	ELISA, WB, IHC
CWF19L1 Polyclonal Antibody	A58694	Human	ELISA, WB, IHC
cwlM Polyclonal Antibody	A64374	Bacillus licheniformis	ELISA
CX3CL1 Polyclonal Antibody	A55627	Human	ELISA, IHC, IF
CXADR Polyclonal Antibody	A58698	Human	ELISA, IHC
Cxcl1 Polyclonal Antibody	A51950	Mouse	ELISA
CXCL1 Polyclonal Antibody	A51978	Human	ELISA
CXCL10 Polyclonal Antibody	A52317	Human	ELISA
CXCL10 Polyclonal Antibody	A56210	Macaca mulatta	ELISA
Cxcl11 Polyclonal Antibody	A54097	Mouse, Human	ELISA, IHC
CXCL13 Polyclonal Antibody	A53209	Human	ELISA, IHC
Cxcl14 Polyclonal Antibody	A53101	Mouse	ELISA
Cxcl16 Polyclonal Antibody	A54204	Mouse	ELISA
Cxcl3 Polyclonal Antibody	A53097	Mouse	ELISA
CXCL3 Polyclonal Antibody	A64116	Human	ELISA, IHC
CXCL5 Polyclonal Antibody	A51689	Human	ELISA, WB
Cxcl5 Polyclonal Antibody	A54093	Mouse, Human	ELISA, WB
CXCL6 Polyclonal Antibody	A58702	Human	ELISA, IHC
CXCL8 Polyclonal Antibody	A51853	Human	ELISA, WB, IHC
CXCL8 Polyclonal Antibody	A52034	Gallus gallus	ELISA
CXCL8 Polyclonal Antibody	A55395	Dog	ELISA
Cxcl9 Polyclonal Antibody	A53125	Mouse	ELISA

Product Name	Cat. #	Species Reactivity	Applications
CXCR2 Polyclonal Antibody	A62782	Human	ELISA, IHC, IF
CXCR4 Polyclonal Antibody	A50165	Human, Mouse	ELISA, WB, IHC
CXorf21 Polyclonal Antibody	A67386	Human	ELISA, IHC
CXorf40A Polyclonal Antibody	A66546	Human	ELISA
CXorf40B Polyclonal Antibody	A66686	Human	ELISA
CXorf56 Polyclonal Antibody	A67714	Human	ELISA, IHC
CXorf58 Polyclonal Antibody	A67114	Human	ELISA
CXXC5 Polyclonal Antibody	A65418	Human	ELISA, IHC
CYB5A Polyclonal Antibody	A51913	Human	ELISA, IF
CYB5B Polyclonal Antibody	A62362	Human, Mouse	ELISA, WB, IHC, IF
CYB5R1 Polyclonal Antibody	A58706	Human	ELISA, IHC
CYB5R2 Polyclonal Antibody	A65362	Human, Mouse	ELISA, WB, IHC
CYB5R3 Polyclonal Antibody	A51645	Human	ELISA, WB, IHC, IF
CYBA Polyclonal Antibody	A68507	Human	ELISA, IF
CYBB Polyclonal Antibody	A50169	Human	ELISA, IHC, IF
CYC1 Polyclonal Antibody	A53893	Human	ELISA, WB, IHC, IF
CYCB1-1 Polyclonal Antibody	A64362	Arabidopsis thaliana	ELISA, WB
CYCS Polyclonal Antibody	A53400	Human	ELISA, IHC
CYGB Polyclonal Antibody	A50173	Human	ELISA, IHC, IF
CYLC2 Polyclonal Antibody	A58710	Human	ELISA, WB
CYP11A1 Polyclonal Antibody	A55579	Human, Rat	ELISA, WB
cyp125 Polyclonal Antibody	A57441	Mycobacterium tuberculosis	ELISA
CYP17A1 Polyclonal Antibody	A54704	Human	ELISA, WB, IHC, IF
CYP17A1 Polyclonal Antibody	A55842	Human	ELISA, WB, IHC, IF
CYP2E1 Polyclonal Antibody	A50177	Human, Mouse	ELISA, WB, IHC
CYP2U1 Polyclonal Antibody	A65794	Human	ELISA, IF
Cyp3a13 Polyclonal Antibody	A65310	Mouse	ELISA
CYP3A7 Polyclonal Antibody	A62366	Human	ELISA, IHC, IF
CYP4F11 Polyclonal Antibody	A51305	Human, Mouse	ELISA, WB, IF
CYP4X1 Polyclonal Antibody	A66842	Human	ELISA, IHC, IF
CYP8B1 Polyclonal Antibody	A68511	Human, Rat	ELISA, WB, IHC, IF
CYS4 Polyclonal Antibody	A56719	Saccharomyces cerevisiae	ELISA
CYSLTR1 Polyclonal Antibody	A62370	Human, Mouse	ELISA, WB, IHC, IF
CYTL1 Polyclonal Antibody	A58714	Human	ELISA, IHC
Cytotoxin 4 Polyclonal Antibody	A64430	Naja atra	ELISA
D2HGDH Polyclonal Antibody	A58718	Human, Mouse	ELISA, WB
d2hgdh Polyclonal Antibody	A64546	Zebrafish	ELISA
Da-3 Polyclonal Antibody	A68767	Trifolium repens	ELISA
DAD1 Polyclonal Antibody	A51601	Human	ELISA, WB, IHC, IF
DALRD3 Polyclonal Antibody	A65122	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
DAND5 Polyclonal Antibody	A56955	Human	ELISA, IHC
DAO Polyclonal Antibody	A50181	Human	ELISA, IHC
DAOA Polyclonal Antibody	A53792	Human, Mouse	ELISA, WB
DAP Polyclonal Antibody	A52804	Human	ELISA, WB
DAPK2 Polyclonal Antibody	A50185	Human, Mouse	ELISA, WB, IHC, IF
DAPL1 Polyclonal Antibody	A62374	Human	ELISA, IHC
DARS Polyclonal Antibody	A62378	Human	ELISA
DAZAP1 Polyclonal Antibody	A66290	Human	ELISA, WB, IF
DBF4B Polyclonal Antibody	A58722	Human	ELISA, WB, IHC
DBI Polyclonal Antibody	A55166	Human	ELISA, IHC
DBNDD2 Polyclonal Antibody	A56823	Human	ELISA, IHC
DCAF4 Polyclonal Antibody	A51189	Human	ELISA, WB
DCAKD Polyclonal Antibody	A58726	Human	ELISA
DCD Polyclonal Antibody	A53288	Human	ELISA
DCN Polyclonal Antibody	A52807	Human, Mouse	ELISA, WB, IHC
Dcn Polyclonal Antibody	A55862	Mouse	ELISA
DCN Polyclonal Antibody	A62386	Bovine	ELISA, WB
DCP2 Polyclonal Antibody	A55910	Human	ELISA, IHC
DCTN2 Polyclonal Antibody	A58730	Human, Mouse	ELISA, WB, IHC
DCTN3 Polyclonal Antibody	A57037	Human	ELISA, IHC
DCTN4 Polyclonal Antibody	A58734	Human, Mouse	ELISA, WB
DCTN5 Polyclonal Antibody	A67442	Human	ELISA, IHC
DCUN1D1 Polyclonal Antibody	A56455	Human	ELISA, IHC
DCUN1D2 Polyclonal Antibody	A65586	Human	ELISA
DCX Polyclonal Antibody	A58738	Human, Mouse	ELISA, WB, IHC, IF
DCX Polyclonal Antibody	A62390	Human	ELISA, IHC, IF
dcyD Polyclonal Antibody	A57669	Escherichia coli	ELISA, WB
DDAH2 Polyclonal Antibody	A62394	Human, Mouse	ELISA, WB, IHC
DDHD2 Polyclonal Antibody	A58742	Human	ELISA, WB, IHC
DDI2 Polyclonal Antibody	A65330	Human	ELISA, WB, IHC
DDIT3 Polyclonal Antibody	A62398	Human	ELISA, IHC, IF
Ddt Polyclonal Antibody	A57381	Mouse	ELISA
DDX17 Polyclonal Antibody	A69103	Human	ELISA, IF
DDX19A Polyclonal Antibody	A55303	Human, Mouse	ELISA, WB, IHC
DDX27 Polyclonal Antibody	A56006	Human	ELISA
DDX28 Polyclonal Antibody	A69235	Human	ELISA, WB
DDX31 Polyclonal Antibody	A58746	Human	ELISA, IHC
DDX39B Polyclonal Antibody	A53845	Human	ELISA, WB, IHC
DDX3X Polyclonal Antibody	A50189	Human, Mouse, Rat	ELISA, WB, IHC, IF
DDX42 Polyclonal Antibody	A58750	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
DDX50 Polyclonal Antibody	A58754	Human	ELISA, WB, IHC
DDX55 Polyclonal Antibody	A64168	Human	ELISA, IHC
DDX58 Polyclonal Antibody	A54640	Human	ELISA
DDX59 Polyclonal Antibody	A65234	Human	ELISA, IHC
def Polyclonal Antibody	A57013	Escherichia coli	ELISA, WB
DEF8 Polyclonal Antibody	A65722	Human	ELISA
Defa5 Polyclonal Antibody	A62402	Mouse	ELISA
DEFB1 Polyclonal Antibody	A52788	Human	ELISA, IHC, IF
DEFB104A Polyclonal Antibody	A66658	Human	ELISA, IHC
DEFB126 Polyclonal Antibody	A57465	Pongo pygmaeus	ELISA
DEFB129 Polyclonal Antibody	A67250	Human	ELISA, WB, IHC
Defb14 Polyclonal Antibody	A55000	Mouse, Rat	ELISA, WB
Defensin D2 Polyclonal Antibody	A64212	Spinacia oleracea	ELISA
DEGP12 Polyclonal Antibody	A51309	Arabidopsis thaliana	ELISA, WB
Delta-Ctenitoxin-Pn1a Polyclonal Antibody	A64406	Phoneutria nigriventer	ELISA
DENND1A Polyclonal Antibody	A58758	Human	ELISA, IHC
DENND2C Polyclonal Antibody	A65246	Human	ELISA, WB
DENND2D Polyclonal Antibody	A58762	Human	ELISA, WB, IHC
DENND5A Polyclonal Antibody	A65602	Human	ELISA, IHC
DENR Polyclonal Antibody	A56022	Human	ELISA
deoD Polyclonal Antibody	A53072	Escherichia coli	ELISA, WB
DEXI Polyclonal Antibody	A58766	Human	ELISA, IHC
DFNA5 Polyclonal Antibody	A55235	Human	ELISA, IHC, IF
DGCR6L Polyclonal Antibody	A57105	Human, Mouse	ELISA, WB
DGKG Polyclonal Antibody	A62406	Human, Mouse	ELISA, WB
DGUOK Polyclonal Antibody	A57313	Human	ELISA, IHC
DHFR Polyclonal Antibody	A50193	Human	ELISA, IHC, IF
DHRS12 Polyclonal Antibody	A58770	Human	ELISA, WB, IHC, IF
DHRS13 Polyclonal Antibody	A58774	Human	ELISA, WB, IHC
DHRS3 Polyclonal Antibody	A62410	Human, Mouse	ELISA, WB, IHC, IF
DHRS7 Polyclonal Antibody	A56847	Human	ELISA, WB, IHC
DHRS7B Polyclonal Antibody	A58778	Human	ELISA
DHRS9 Polyclonal Antibody	A53477	Human	ELISA, IHC, IF
DHRSX Polyclonal Antibody	A58782	Human	ELISA, IHC
DHX15 Polyclonal Antibody	A68515	Human	ELISA, WB
DHX30 Polyclonal Antibody	A58786	Human	ELISA, IHC
DHX32 Polyclonal Antibody	A65494	Human, Mouse	ELISA, WB, IHC
DHX34 Polyclonal Antibody	A69895	Human	ELISA, IHC, IF
DHX35 Polyclonal Antibody	A67950	Human, Mouse	ELISA, WB, IHC
DHX9 Polyclonal Antibody	A56274	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
DIMT1 Polyclonal Antibody	A56992	Human, Mouse	ELISA, WB, IHC
DIO1 Polyclonal Antibody	A64386	Human	ELISA
DIO2 Polyclonal Antibody	A70027	Human	ELISA, IHC, IF
Dio3 Polyclonal Antibody	A55826	Rat	ELISA
DIP2A Polyclonal Antibody	A58790	Human	ELISA, IHC
DIRC1 Polyclonal Antibody	A66814	Human	ELISA, IHC
DIS3L Polyclonal Antibody	A58794	Human	ELISA, WB, IHC
DIS3L2 Polyclonal Antibody	A58798	Human	ELISA, IHC
Disintegrin EC3B Polyclonal Antibody	A54974	Echis carinatus	ELISA
DKK1 Polyclonal Antibody	A64088	Human	ELISA, IHC, IF
DKK2 Polyclonal Antibody	A56671	Human	ELISA
DKK3 Polyclonal Antibody	A58802	Human	ELISA
DKKL1 Polyclonal Antibody	A68134	Human	ELISA, IHC
DLEU1 Polyclonal Antibody	A58806	Human	ELISA, IHC, IF
DLK1 Polyclonal Antibody	A54519	Human	ELISA, WB, IHC, IF
Dlk2 Polyclonal Antibody	A66110	Mouse	ELISA, WB
DLL3 Polyclonal Antibody	A55782	Human	ELISA
DLST Polyclonal Antibody	A69175	Human, Rat	ELISA, WB, IHC, IF
DLX1 Polyclonal Antibody	A68519	Human	ELISA, IF
DMGDH Polyclonal Antibody	A69267	Human, Mouse	ELISA, WB, IF
DMKN Polyclonal Antibody	A58810	Human, Mouse	ELISA, WB
DMRT3 Polyclonal Antibody	A70315	Human	ELISA, IHC
DMRTA1 Polyclonal Antibody	A65294	Human	ELISA, IHC
DNA2 Polyclonal Antibody	A62414	Human	ELISA, IHC
DNAAF5 Polyclonal Antibody	A58814	Human	ELISA
DNAH3 Polyclonal Antibody	A56887	Human	ELISA, IHC
DNAJB1 Polyclonal Antibody	A51817	Human, Mouse	ELISA, WB, IHC
DNAJB14 Polyclonal Antibody	A66234	Human	ELISA, IHC, IF
DNAJB2 Polyclonal Antibody	A57017	Human	ELISA, WB, IHC
DNAJB4 Polyclonal Antibody	A50668	Human, Mouse	ELISA, WB, IHC
DNAJB5 Polyclonal Antibody	A50197	Human	ELISA, WB, IHC
DNAJB8 Polyclonal Antibody	A58818	Human, Mouse	ELISA, WB, IHC
DNAJC17 Polyclonal Antibody	A58822	Human	ELISA, WB, IHC
DNAJC30 Polyclonal Antibody	A56807	Human	ELISA, WB, IHC
DNAJC6 Polyclonal Antibody	A62418	Human	ELISA, IHC, IF
DNAL4 Polyclonal Antibody	A62422	Human, Mouse	ELISA, WB, IHC, IF
DNALI1 Polyclonal Antibody	A62426	Human, Mouse	ELISA, WB, IHC, IF
DNASE1L3 Polyclonal Antibody	A64790	Human	ELISA, IHC
DNM1 Polyclonal Antibody	A50201	Human	ELISA, IHC, IF
DNM1L Polyclonal Antibody	A62430	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
DNMT3A Polyclonal Antibody	A52569	Human, Mouse	ELISA, IHC, IF
DNPEP Polyclonal Antibody	A69463	Human	ELISA, IHC, IF
DNTT Polyclonal Antibody	A58826	Human	ELISA, IHC, IF
DNTTIP1 Polyclonal Antibody	A54220	Human	ELISA, IHC
DOCK10 Polyclonal Antibody	A58830	Human	ELISA, IHC
DOCK5 Polyclonal Antibody	A67574	Human	ELISA
DODA Polyclonal Antibody	A69955	Portulaca grandiflora	ELISA
DOK1 Polyclonal Antibody	A55198	Human	ELISA, IHC, IF
DOK5 Polyclonal Antibody	A54752	Human, Mouse	ELISA, WB, IHC
DOK6 Polyclonal Antibody	A58834	Human	ELISA
DPH1 Polyclonal Antibody	A51301	Human	ELISA, IHC
DPH2 Polyclonal Antibody	A51225	Human, Mouse	ELISA, WB, IHC
DPH3 Polyclonal Antibody	A58838	Human	ELISA
DPH6 Polyclonal Antibody	A65542	Human	ELISA, IHC, IF
DPP6 Polyclonal Antibody	A54660	Human	ELISA
DPPA3 Polyclonal Antibody	A68959	Human	ELISA, IHC
DPPA4 Polyclonal Antibody	A58842	Human	ELISA, IHC
DPPA5 Polyclonal Antibody	A69467	Human, Mouse	ELISA, WB, IHC, IF
DPY19L1 Polyclonal Antibody	A69811	Human	ELISA, IF
DPY19L3 Polyclonal Antibody	A69947	Human	ELISA, WB, IF
DPY30 Polyclonal Antibody	A58846	Human	ELISA, IHC
DRAP1 Polyclonal Antibody	A51849	Human	ELISA
DRC7 Polyclonal Antibody	A58850	Human	ELISA, IHC
DRD5 Polyclonal Antibody	A50205	Human, Mouse	ELISA, WB, IHC, IF
DRG2 Polyclonal Antibody	A58854	Human	ELISA
dsbA Polyclonal Antibody	A50881	Pseudomonas syringae pv. tomato	ELISA
dsbA Polyclonal Antibody	A57229	Pseudomonas syringae pv. tomato	ELISA
DSC3 Polyclonal Antibody	A69847	Human	ELISA, IHC, IF
DSCAM Polyclonal Antibody	A69471	Human	ELISA, IHC
DSCC1 Polyclonal Antibody	A67918	Human	ELISA, WB
DSCR4 Polyclonal Antibody	A62434	Human	ELISA
DSCR9 Polyclonal Antibody	A62438	Human	ELISA, IHC
dsdA Polyclonal Antibody	A56394	Salmonella heidelberg	ELISA
dsdA Polyclonal Antibody	A56439	Salmonella heidelberg	ELISA
DSE Polyclonal Antibody	A58858	Human	ELISA, IHC, IF
DSG3 Polyclonal Antibody	A54910	Human	ELISA, IHC
Dsg3 Polyclonal Antibody	A58862	Mouse	ELISA
DSP Polyclonal Antibody	A62442	Human	ELISA
DSTN Polyclonal Antibody	A58866	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
DTNBP1 Polyclonal Antibody	A69059	Human	ELISA, IHC, IF
DTWD1 Polyclonal Antibody	A53388	Human, Mouse	ELISA, IHC, IF
DUS3L Polyclonal Antibody	A58870	Human	ELISA, IHC
DUS4L Polyclonal Antibody	A62446	Human	ELISA, IHC
DUSP13 Polyclonal Antibody	A55339	Human	ELISA, WB, IHC
DUSP19 Polyclonal Antibody	A67062	Human	ELISA, IHC, IF
DUSP21 Polyclonal Antibody	A67570	Human	ELISA, IHC
DUSP26 Polyclonal Antibody	A51241	Human, Mouse	ELISA, WB, IHC
DYDC1 Polyclonal Antibody	A66262	Human	ELISA, WB
DYNLL1 Polyclonal Antibody	A53885	Human	ELISA, WB, IHC, IF
DYNLL2 Polyclonal Antibody	A58874	Human	ELISA
DYNLRB1 Polyclonal Antibody	A51253	Human	ELISA, IHC
DYRK3 Polyclonal Antibody	A62450	Human	ELISA
DYRK4 Polyclonal Antibody	A58878	Human	ELISA, IHC, IF
E2 Polyclonal Antibody	A55218	Human papillomavirus type 18	ELISA
E2F2 Polyclonal Antibody	A50937	Human	ELISA, WB, IHC
E6 Polyclonal Antibody	A55631	Human papillomavirus type 16	ELISA
E6 Polyclonal Antibody	A64350	Human papillomavirus type 52	ELISA
E7 Polyclonal Antibody	A55263	Human papillomavirus type 18	ELISA
EAPP Polyclonal Antibody	A58882	Human, Mouse	ELISA, WB, IHC
EBF3 Polyclonal Antibody	A69251	Human	ELISA, WB
EBI3 Polyclonal Antibody	A62454	Human	ELISA, WB, IHC
EBP Polyclonal Antibody	A50941	Human	ELISA, WB, IHC, IF
ECD Polyclonal Antibody	A58886	Human	ELISA, WB
ECHDC1 Polyclonal Antibody	A67738	Human	ELISA
ECM1 Polyclonal Antibody	A69475	Human	ELISA, WB, IHC
ecoRVM Polyclonal Antibody	A57405	Escherichia coli	ELISA
EDC4 Polyclonal Antibody	A68951	Human	ELISA, IHC, IF
EDEM2 Polyclonal Antibody	A58890	Human	ELISA, IHC
EEF1D Polyclonal Antibody	A54224	Human	ELISA, WB
EEF1E1 Polyclonal Antibody	A55990	Human	ELISA, WB, IHC, IF
EEF2 Polyclonal Antibody	A53432	Human	ELISA, IHC, IF
EEF2K Polyclonal Antibody	A62458	Human	ELISA, IHC, IF
EEF2KMT Polyclonal Antibody	A58894	Human	ELISA, WB, IHC
EFCAB1 Polyclonal Antibody	A67202	Human	ELISA
EFCAB12 Polyclonal Antibody	A65382	Human	ELISA, IHC
EFCAB14 Polyclonal Antibody	A58898	Human	ELISA, WB, IHC
EFHB Polyclonal Antibody	A58902	Human	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
Efna1 Polyclonal Antibody	A53384	Rat	ELISA
Efna5 Polyclonal Antibody	A53141	Rat	ELISA
Efnb3 Polyclonal Antibody	A56631	Mouse, Human	ELISA, WB
EFR3B Polyclonal Antibody	A68206	Human	ELISA, IHC
EGF Polyclonal Antibody	A52110	Human	ELISA
EGFL6 Polyclonal Antibody	A58906	Human	ELISA, WB, IHC
EGFLAM Polyclonal Antibody	A65242	Human, Mouse	ELISA, WB
EGFR Polyclonal Antibody	A52964	Human	ELISA, IHC
EGLN2 Polyclonal Antibody	A54934	Human	ELISA, IHC, IF
Egln3 Polyclonal Antibody	A52097	Mouse	ELISA
EGR1 Polyclonal Antibody	A54559	Human	ELISA
EGR3 Polyclonal Antibody	A70359	Human	ELISA, WB
egsA Polyclonal Antibody	A64566	Geobacillus stearothermophilus	ELISA
EHBP1 Polyclonal Antibody	A66510	Human	ELISA
EHHADH Polyclonal Antibody	A56747	Human, Mouse	ELISA, WB, IHC
EIF1AD Polyclonal Antibody	A66958	Human	ELISA
EIF3C Polyclonal Antibody	A67130	Human, Mouse	ELISA, WB, IHC, IF
EIF3D Polyclonal Antibody	A62554	Human, Mouse, Rat	ELISA, WB, IHC
EIF3G Polyclonal Antibody	A50620	Human	ELISA, WB
EIF3I Polyclonal Antibody	A51665	Human	ELISA, WB
EIF4A2 Polyclonal Antibody	A52667	Human	ELISA, WB, IHC
EIF4A2 Polyclonal Antibody	A53857	Human	ELISA, WB
EIF4EBP2 Polyclonal Antibody	A53121	Human	ELISA, IHC
EIF4EBP2 Polyclonal Antibody	A53555	Human, Mouse	ELISA, WB, IHC
EIF4EBP3 Polyclonal Antibody	A50209	Human	ELISA, IHC
EIF5 Polyclonal Antibody	A56082	Human	ELISA, IHC, IF
EIF5A2 Polyclonal Antibody	A67474	Human	ELISA, IHC, IF
EIF6 Polyclonal Antibody	A62462	Human	ELISA, IHC, IF
ELAC1 Polyclonal Antibody	A58910	Human	ELISA, IHC
ELANE Polyclonal Antibody	A56090	Human	ELISA, IHC, IF
ELAVL1 Polyclonal Antibody	A58914	Human	ELISA, WB
ELAVL2 Polyclonal Antibody	A53190	Human, Mouse	ELISA, WB, IHC, IF
ELAVL4 Polyclonal Antibody	A58918	Human	ELISA, WB, IHC, IF
ELMOD3 Polyclonal Antibody	A70091	Human	ELISA, WB, IHC, IF
ELOVL1 Polyclonal Antibody	A58922	Human	ELISA, IHC, IF
ELOVL2 Polyclonal Antibody	A67990	Human	ELISA, IHC, IF
ELOVL5 Polyclonal Antibody	A67758	Human	ELISA, IF
ELP3 Polyclonal Antibody	A69279	Human, Rat	ELISA, WB, IF
eltB Polyclonal Antibody	A55355	Escherichia coli	ELISA
EMC6 Polyclonal Antibody	A58926	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
EMC8 Polyclonal Antibody	A58930	Human	ELISA, WB, IHC
EMC9 Polyclonal Antibody	A57257	Human	ELISA
EME1 Polyclonal Antibody	A51209	Human	ELISA, IHC
EMILIN1 Polyclonal Antibody	A58934	Human	ELISA, IHC
EML2 Polyclonal Antibody	A56010	Human	ELISA, IHC, IF
EMP1 Polyclonal Antibody	A62466	Human	ELISA, IHC, IF
EMP3 Polyclonal Antibody	A62470	Human, Mouse	ELISA, WB, IF
ENC1 Polyclonal Antibody	A58938	Human, Mouse	ELISA, WB
ENKD1 Polyclonal Antibody	A58942	Human, Mouse	ELISA, WB, IHC
ENO1 Polyclonal Antibody	A52700	Human, Mouse	ELISA, WB, IHC, IF
Eno1 Polyclonal Antibody	A52768	Human, Mouse	ELISA, WB
ENO2 Polyclonal Antibody	A51942	Human, Mouse	ELISA, WB, IHC
ENO2 Polyclonal Antibody	A54248	Human	ELISA
ENO3 Polyclonal Antibody	A52286	Human	ELISA, IHC, IF
ENOSF1 Polyclonal Antibody	A69939	Human	ELISA, IHC
ENOX1 Polyclonal Antibody	A62474	Human	ELISA, WB, IHC
ENPEP Polyclonal Antibody	A52154	Human	ELISA, IHC
ENPP6 Polyclonal Antibody	A65390	Human	ELISA, IHC, IF
ENSA Polyclonal Antibody	A57615	Human	ELISA, IHC, IF
entC1 Polyclonal Antibody	A50849	Staphylococcus aureus	ELISA
entC2 Polyclonal Antibody	A64370	Staphylococcus aureus	ELISA
entC3 Polyclonal Antibody	A50825	Staphylococcus aureus	ELISA
entD Polyclonal Antibody	A50776	Staphylococcus aureus	ELISA
entE Polyclonal Antibody	A50744	Staphylococcus aureus	ELISA
entG Polyclonal Antibody	A50841	Staphylococcus aureus	ELISA
entG Polyclonal Antibody	A64478	Staphylococcus aureus	ELISA
entH Polyclonal Antibody	A64482	Staphylococcus aureus	ELISA
ENTPD3 Polyclonal Antibody	A58946	Human	ELISA, IHC, IF
ENTPD7 Polyclonal Antibody	A62478	Human, Mouse	ELISA, WB, IF
ENTPD8 Polyclonal Antibody	A62482	Human, Mouse	ELISA, WB, IHC, IF
ENY2 Polyclonal Antibody	A58950	Human	ELISA
EPB2 Polyclonal Antibody	A64310	Hordeum vulgare	ELISA
EPCAM Polyclonal Antibody	A68759	Human, Rat	ELISA, WB, IHC, IF
EPHA7 Polyclonal Antibody	A64722	Human	ELISA, IHC
EPHX1 Polyclonal Antibody	A56373	Human	ELISA, WB, IHC
EPHX2 Polyclonal Antibody	A50213	Human	ELISA, WB, IHC
EPM2AIP1 Polyclonal Antibody	A65634	Human	ELISA, IHC
EPO Polyclonal Antibody	A53364	Human	ELISA, IHC
EPO Polyclonal Antibody	A62486	Bovine	ELISA, WB
EPO Polyclonal Antibody	A62487	Macaca fascicularis	ELISA

Product Name	Cat. #	Species Reactivity	Applications
EPO Polyclonal Antibody	A64092	Human, Mouse	ELISA, WB, IHC, IF
EPPK1 Polyclonal Antibody	A54372	Human	ELISA
EPRS Polyclonal Antibody	A69479	Human	ELISA, WB, IHC, IF
EPS8L2 Polyclonal Antibody	A67958	Human	ELISA, IHC
EPX Polyclonal Antibody	A69483	Human	ELISA, IHC
ERAS Polyclonal Antibody	A52309	Human	ELISA, WB
ERBB2 Polyclonal Antibody	A52999	Human	ELISA, IHC
ERCC6L2 Polyclonal Antibody	A62190	Human	ELISA, IHC
ERG Polyclonal Antibody	A62498	Human, Rat	ELISA, IP
ERGIC1 Polyclonal Antibody	A51109	Human	ELISA, WB, IHC, IF
ERGIC3 Polyclonal Antibody	A55615	Human	ELISA, WB, IHC, IF
ERH Polyclonal Antibody	A51685	Human	ELISA, IHC
ERLIN1 Polyclonal Antibody	A58954	Human, Mouse	ELISA, WB, IHC
ERO1B Polyclonal Antibody	A65938	Human, Mouse	ELISA, WB, IHC
ERP27 Polyclonal Antibody	A58958	Human	ELISA, IHC
ERP29 Polyclonal Antibody	A51705	Human	ELISA, WB, IHC
ERP29 Polyclonal Antibody	A51725	Human	ELISA, WB, IHC
ERVK-6 Polyclonal Antibody	A65250	Human	ELISA
ESAM Polyclonal Antibody	A55335	Human	ELISA, IHC
ESD Polyclonal Antibody	A62502	Human	ELISA, IF
ESM1 Polyclonal Antibody	A58962	Human	ELISA
ESR1 Polyclonal Antibody	A55755	Human	ELISA, IHC, IF
ESR1 Polyclonal Antibody	A68315	Human, Rat	ELISA, WB, IHC, ChIP
ESR2 Polyclonal Antibody	A53738	Human, Mouse	ELISA, WB, IHC
ESRG Polyclonal Antibody	A58966	Human	ELISA, IHC
ESRRB Polyclonal Antibody	A69487	Human	ELISA, WB
ESRRG Polyclonal Antibody	A62506	Human	ELISA
esxB Polyclonal Antibody	A57738	Mycobacterium tuberculosis	ELISA
ESYT2 Polyclonal Antibody	A68523	Human	ELISA, IF
ETFDH Polyclonal Antibody	A55539	Pig	ELISA
ETFDH Polyclonal Antibody	A58970	Human, Mouse	ELISA, WB, IHC
ETHE1 Polyclonal Antibody	A58974	Human, Mouse	ELISA, WB, IHC
ETNK1 Polyclonal Antibody	A58978	Human	ELISA, WB, IHC, IF
ETS1 Polyclonal Antibody	A62510	Human	ELISA
eutC Polyclonal Antibody	A52496	Escherichia coli	ELISA, WB
EVA1C Polyclonal Antibody	A58982	Human	ELISA, WB, IHC
EVI2A Polyclonal Antibody	A62514	Human	ELISA, IHC, IF
EXO1 Polyclonal Antibody	A62518	Human	ELISA, WB, IHC
EXOSC1 Polyclonal Antibody	A68214	Human	ELISA, IHC
EXOSC3 Polyclonal Antibody	A67974	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
EXOSC5 Polyclonal Antibody	A58986	Human	ELISA, WB, IHC
EXOSC8 Polyclonal Antibody	A66990	Human, Mouse	ELISA, WB
EXTL3 Polyclonal Antibody	A51669	Human	ELISA, WB, IHC, IF
EYA1 Polyclonal Antibody	A69151	Human	ELISA, IF
EYA2 Polyclonal Antibody	A55675	Human	ELISA, IHC
EYA4 Polyclonal Antibody	A69491	Human	ELISA, IHC, IF
EZR Polyclonal Antibody	A55547	Human	ELISA
F11 Polyclonal Antibody	A62522	Human	ELISA, IHC
F2 Polyclonal Antibody	A58990	Human	ELISA, WB, IF
F2R Polyclonal Antibody	A62526	Human	ELISA, IHC, IF
F3 Polyclonal Antibody	A50217	Human, Mouse	ELISA, WB, IHC, IF
F7 Polyclonal Antibody	A62530	Human	ELISA, IHC
F8 Polyclonal Antibody	A62534	Human	ELISA, IHC, IF
fabD Polyclonal Antibody	A50961	Staphylococcus aureus	ELISA
fabF Polyclonal Antibody	A64502	Escherichia coli	ELISA
fabF Polyclonal Antibody	A65102	Staphylococcus aureus	ELISA
fabI Polyclonal Antibody	A65498	Staphylococcus aureus, Mouse	ELISA, WB
FABP1 Polyclonal Antibody	A54672	Human	ELISA, WB, IHC
FABP3 Polyclonal Antibody	A52893	Human, Mouse, Rat	ELISA, WB, IHC
FABP3 Polyclonal Antibody	A54216	Human	ELISA
FABP3 Polyclonal Antibody	A54252	Human	ELISA
FAH Polyclonal Antibody	A62538	Human, Mouse	ELISA, WB, IHC, IF
FAHD2A Polyclonal Antibody	A66294	Human	ELISA, WB, IHC
FAIM Polyclonal Antibody	A58994	Human, Rat	ELISA, WB, IHC
FAM102B Polyclonal Antibody	A65350	Human	ELISA, IHC, IF
FAM105A Polyclonal Antibody	A58998	Human	ELISA, WB, IHC
FAM107A Polyclonal Antibody	A62542	Human	ELISA, IHC
FAM107B Polyclonal Antibody	A59002	Human, Mouse	ELISA, WB
FAM109B Polyclonal Antibody	A65554	Human	ELISA
FAM110B Polyclonal Antibody	A53567	Human, Mouse	ELISA, WB, IHC
FAM114A2 Polyclonal Antibody	A55731	Human, Mouse	ELISA, WB, IHC
FAM117B Polyclonal Antibody	A65578	Human	ELISA, IHC
FAM118A Polyclonal Antibody	A59006	Human	ELISA, WB, IHC, IF
FAM118B Polyclonal Antibody	A59010	Human	ELISA, WB, IHC
FAM122B Polyclonal Antibody	A65782	Human, Mouse	ELISA, WB, IHC
FAM124A Polyclonal Antibody	A65958	Human	ELISA, IHC
FAM124B Polyclonal Antibody	A59014	Human	ELISA, WB, IHC
FAM126B Polyclonal Antibody	A59018	Human	ELISA, WB, IHC
FAM127A Polyclonal Antibody	A62546	Human	ELISA, WB, IHC
FAM127B Polyclonal Antibody	A59022	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
FAM127C Polyclonal Antibody	A64806	Human	ELISA, IHC
FAM131A Polyclonal Antibody	A59026	Human	ELISA, IHC
FAM131B Polyclonal Antibody	A65930	Human	ELISA
FAM135B Polyclonal Antibody	A65022	Human	ELISA, IHC
FAM136A Polyclonal Antibody	A59030	Human	ELISA, WB, IHC
FAM149B1 Polyclonal Antibody	A59034	Human, Mouse	ELISA, WB, IHC
FAM151B Polyclonal Antibody	A65398	Human, Mouse	ELISA, WB, IHC, IF
FAM153C Polyclonal Antibody	A65006	Human	ELISA, IHC
FAM156A Polyclonal Antibody	A59038	Human, Mouse	ELISA, WB, IHC, IF
FAM160A2 Polyclonal Antibody	A66374	Human	ELISA, IHC
FAM160B1 Polyclonal Antibody	A55942	Human	ELISA, IHC
FAM161A Polyclonal Antibody	A64930	Human, Mouse	ELISA, WB, IHC
FAM163A Polyclonal Antibody	A59042	Human	ELISA, IHC
FAM166A Polyclonal Antibody	A65562	Human	ELISA, IHC
FAM167A Polyclonal Antibody	A66462	Human	ELISA, IHC
FAM168B Polyclonal Antibody	A62550	Human	ELISA, IHC, IF
FAM171B Polyclonal Antibody	A65610	Human	ELISA, IHC, IF
FAM174A Polyclonal Antibody	A59046	Human	ELISA, IHC
FAM182A Polyclonal Antibody	A65274	Human	ELISA, IF
FAM187B Polyclonal Antibody	A64834	Human	ELISA, IHC
FAM19A1 Polyclonal Antibody	A65830	Human	ELISA
FAM19A4 Polyclonal Antibody	A59050	Human	ELISA
FAM20C Polyclonal Antibody	A70015	Human	ELISA, WB, IHC
FAM222B Polyclonal Antibody	A66978	Human, Mouse	ELISA, WB, IHC
FAM234B Polyclonal Antibody	A62826	Human	ELISA, WB, IHC, IF
FAM32A Polyclonal Antibody	A59054	Human	ELISA, IHC
FAM3B Polyclonal Antibody	A62562	Human, Mouse, Rat	ELISA, WB, IHC
FAM3C Polyclonal Antibody	A67078	Human	ELISA, WB
FAM43A Polyclonal Antibody	A67034	Human, Mouse	ELISA, WB, IHC
FAM45A Polyclonal Antibody	A66650	Human	ELISA
FAM46B Polyclonal Antibody	A59058	Human	ELISA, WB
FAM49B Polyclonal Antibody	A57265	Human	ELISA
FAM53B Polyclonal Antibody	A59062	Human	ELISA, WB, IHC
FAM60A Polyclonal Antibody	A51357	Human	ELISA, IHC
FAM63A Polyclonal Antibody	A66742	Human, Mouse	ELISA, WB
FAM69A Polyclonal Antibody	A65326	Human, Mouse	ELISA, WB
FAM71A Polyclonal Antibody	A59066	Human	ELISA, IHC, IF
FAM71C Polyclonal Antibody	A59070	Human	ELISA, WB
FAM71E2 Polyclonal Antibody	A59074	Human	ELISA
FAM71F1 Polyclonal Antibody	A67022	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
FAM71F2 Polyclonal Antibody	A65514	Human	ELISA, IHC
FAM76B Polyclonal Antibody	A59078	Human	ELISA
FAM78A Polyclonal Antibody	A65154	Human	ELISA, WB, IHC
FAM84A Polyclonal Antibody	A70103	Human	ELISA, IHC
FAM89B Polyclonal Antibody	A67042	Human	ELISA
FAM92A1 Polyclonal Antibody	A59082	Human, Mouse	ELISA, WB, IHC, IF
FAM96A Polyclonal Antibody	A59086	Human, Mouse	ELISA, WB
FANCC Polyclonal Antibody	A62566	Human, Mouse	ELISA, WB, IHC, IF
FANCD2 Polyclonal Antibody	A67350	Human	ELISA, IHC
FANCL Polyclonal Antibody	A69263	Human	ELISA, IF
FANK1 Polyclonal Antibody	A56038	Human	ELISA, IHC
FAR2 Polyclonal Antibody	A66314	Human	ELISA, IHC, IF
FARP1 Polyclonal Antibody	A68226	Human	ELISA, WB, IHC, IF
FARP2 Polyclonal Antibody	A59090	Human	ELISA, IHC, IF
FARSA Polyclonal Antibody	A55531	Human	ELISA, IHC, IF
FAS Polyclonal Antibody	A53081	Human	ELISA, WB, IHC
FASLG Polyclonal Antibody	A55854	Human	ELISA, IHC
FASN Polyclonal Antibody	A50700	Human	ELISA, IHC
FASTKD5 Polyclonal Antibody	A69951	Human	ELISA, IHC
FATE1 Polyclonal Antibody	A66582	Human	ELISA, WB, IF
FBLN2 Polyclonal Antibody	A69495	Human	ELISA, WB, IHC
FBLN7 Polyclonal Antibody	A65082	Human	ELISA, IHC
fbpA Polyclonal Antibody	A57689	Mycobacterium tuberculosis	ELISA, WB
FBXL17 Polyclonal Antibody	A67334	Human	ELISA, IHC
FBXL18 Polyclonal Antibody	A54902	Human	ELISA, IHC, IF
FBXL2 Polyclonal Antibody	A59094	Human	ELISA
FBXL20 Polyclonal Antibody	A70031	Human	ELISA, IHC
FBXO15 Polyclonal Antibody	A59098	Human	ELISA, IHC
FBXO24 Polyclonal Antibody	A62570	Human, Mouse	ELISA, WB
FBXO25 Polyclonal Antibody	A66882	Human, Mouse	ELISA, WB, IHC, IF
FBXO28 Polyclonal Antibody	A67982	Human	ELISA, IHC
FBXO30 Polyclonal Antibody	A59102	Human, Mouse	ELISA, WB, IHC
FBXO40 Polyclonal Antibody	A67774	Human	ELISA, IHC
FBXO44 Polyclonal Antibody	A67498	Human	ELISA
FBXW10 Polyclonal Antibody	A59106	Human	ELISA, IHC
FBXW12 Polyclonal Antibody	A65530	Human	ELISA, IHC
FBXW2 Polyclonal Antibody	A68070	Human, Mouse	ELISA, WB, IHC
FCER1G Polyclonal Antibody	A54646	Human	ELISA, IHC, IF
FCER2 Polyclonal Antibody	A56358	Human	ELISA
Fcgr3 Polyclonal Antibody	A59110	Mouse	ELISA

Product Name	Cat. #	Species Reactivity	Applications
FCGRT Polyclonal Antibody	A53808	Human, Mouse	ELISA, WB, IHC, IF
FCGRT Polyclonal Antibody	A55790	Macaca fascicularis	ELISA
FCHO1 Polyclonal Antibody	A69499	Human	ELISA, IHC, IF
FCHSD2 Polyclonal Antibody	A59114	Human	ELISA, WB, IHC
Fcmr Polyclonal Antibody	A56202	Mouse, Rat	ELISA, WB
FCRL4 Polyclonal Antibody	A59118	Human, Mouse	ELISA, WB
FCRL6 Polyclonal Antibody	A65198	Human	ELISA
FDFT1 Polyclonal Antibody	A62574	Human	ELISA, WB, IHC, IF
FDXR Polyclonal Antibody	A59122	Human, Mouse	ELISA, WB, IHC
FERD3L Polyclonal Antibody	A66934	Human	ELISA, IHC, IF
FERMT1 Polyclonal Antibody	A70263	Human	ELISA, IHC, IF
FERMT2 Polyclonal Antibody	A70155	Human	ELISA, IHC, IF
FFAR1 Polyclonal Antibody	A68527	Human	ELISA, IHC, IF
FFAR2 Polyclonal Antibody	A62578	Human	ELISA, IHC, IF
FFAR3 Polyclonal Antibody	A62582	Human, Mouse	ELISA, WB, IHC, IF
FGA Polyclonal Antibody	A56318	Human	ELISA, IHC
FGD6 Polyclonal Antibody	A65666	Human	ELISA, WB, IHC
Fgf1 Polyclonal Antibody	A52780	Mouse, Human	ELISA, WB, IHC
FGF1 Polyclonal Antibody	A55415	Human	ELISA
FGF13 Polyclonal Antibody	A59126	Human, Mouse	ELISA, WB, IHC
FGF2 Polyclonal Antibody	A50696	Human	ELISA, IHC
FGF2 Polyclonal Antibody	A52198	Mouse, Human	ELISA, WB
FGF2 Polyclonal Antibody	A52313	Human, Mouse	ELISA, WB, IHC
FGF21 Polyclonal Antibody	A52640	Human	ELISA, WB, IHC
Fgf21 Polyclonal Antibody	A62586	Mouse	ELISA
Fgf23 Polyclonal Antibody	A55822	Rat	ELISA, WB
FGF7 Polyclonal Antibody	A56430	Human	ELISA
FGFBP2 Polyclonal Antibody	A51373	Human	ELISA, WB, IHC
FGFR1OP2 Polyclonal Antibody	A62590	Human, Mouse	ELISA, WB
FGFR3 Polyclonal Antibody	A53797	Human	ELISA, WB
FGFRL1 Polyclonal Antibody	A68531	Human	ELISA, IHC, IF
FGL2 Polyclonal Antibody	A50925	Human	ELISA, IHC
FH Polyclonal Antibody	A62594	Human, Mouse	ELISA, WB, IHC, IF
Fhit Polyclonal Antibody	A57213	Mouse	ELISA
FHIT Polyclonal Antibody	A57805	Human	ELISA, WB
FIBIN Polyclonal Antibody	A56170	Human	ELISA
FIBP Polyclonal Antibody	A50225	Human	ELISA, WB, IHC, IF
FICD Polyclonal Antibody	A70283	Human, Rat	ELISA, WB, IHC, IF
FIGF Polyclonal Antibody	A53236	Human	ELISA
FILIP1L Polyclonal Antibody	A59130	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
fim2 Polyclonal Antibody	A54740	Bordetella pertussis	ELISA
FITC Polyclonal Antibody	A57819		ELISA
FKBP11 Polyclonal Antibody	A59134	Human, Mouse	ELISA, WB, IHC, IF
FKBP14 Polyclonal Antibody	A67318	Human, Mouse	ELISA, WB
FKBP1A Polyclonal Antibody	A51541	Human	ELISA, WB, IHC, IF
FKBP3 Polyclonal Antibody	A56475	Human, Mouse	ELISA, WB, IHC, IF
FKBP6 Polyclonal Antibody	A68278	Human	ELISA, IHC, IF
FKBP9P1 Polyclonal Antibody	A59138	Human	ELISA
fkpA Polyclonal Antibody	A64426	Escherichia coli O6:H1	ELISA
FKRP Polyclonal Antibody	A70191	Human	ELISA, IHC
fla Polyclonal Antibody	A57084	Borrelia burgdorferi	ELISA
FLAD1 Polyclonal Antibody	A51185	Human	ELISA, WB, IHC
FLG Polyclonal Antibody	A55499	Human	ELISA, IHC
fliC Polyclonal Antibody	A53224	Escherichia coli	ELISA
FLII Polyclonal Antibody	A68843	Human	ELISA, IF
Flii Polyclonal Antibody	A54874	Mouse	ELISA
FLRT1 Polyclonal Antibody	A51261	Human, Mouse	ELISA, WB, IHC
FLT3LG Polyclonal Antibody	A57137	Human	ELISA, IHC
FMNL1 Polyclonal Antibody	A59142	Human	ELISA, IHC
FMO4 Polyclonal Antibody	A62602	Human	ELISA, IHC, IF
FN1 Polyclonal Antibody	A52821	Human	ELISA
FNDC3B Polyclonal Antibody	A59146	Human	ELISA, IHC, IF
FNDC4 Polyclonal Antibody	A67502	Human	ELISA
FNDC7 Polyclonal Antibody	A65354	Human	ELISA
FNTB Polyclonal Antibody	A62606	Human	ELISA, WB, IHC, IF, IP
FOLH1 Polyclonal Antibody	A54153	Human	ELISA, WB, IHC
FOSL2 Polyclonal Antibody	A55327	Human	ELISA, IHC
FOXL2 Polyclonal Antibody	A69503	Human	ELISA, WB
FOXP1 Polyclonal Antibody	A68303	Human	ELISA, WB, IHC, ChIP
FOXP3 Polyclonal Antibody	A52825	Human, Mouse	ELISA, WB, IHC
FOXR2 Polyclonal Antibody	A65622	Human	ELISA, WB, IHC
FPGT Polyclonal Antibody	A53828	Human	ELISA
FPR1 Polyclonal Antibody	A59150	Human	ELISA, WB, IHC, IF
FPR2 Polyclonal Antibody	A59154	Human	ELISA, WB, IF
fre Polyclonal Antibody	A57902	Escherichia coli	ELISA, WB
FREM2 Polyclonal Antibody	A68931	Human	ELISA, IF
FRS3 Polyclonal Antibody	A62610	Human	ELISA, IHC
FSCB Polyclonal Antibody	A59158	Human, Mouse	ELISA, WB, IHC
FSCN2 Polyclonal Antibody	A62614	Human	ELISA
FSHB Polyclonal Antibody	A64148	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
FSIP1 Polyclonal Antibody	A66766	Human	ELISA, IHC
FTH1 Polyclonal Antibody	A51982	Human	ELISA
FTH1 Polyclonal Antibody	A53909	Human	ELISA
FTL Polyclonal Antibody	A52585	Human	ELISA, IHC, IF
FTL Polyclonal Antibody	A55699	Human	ELISA, IHC
FTL3 Polyclonal Antibody	A68763	Soybean	ELISA
FTSJ1 Polyclonal Antibody	A53610	Human, Mouse	ELISA, WB, IHC
FUCA2 Polyclonal Antibody	A52174	Human	ELISA, WB, IHC
FUK Polyclonal Antibody	A70107	Human	ELISA, IHC, IF
FUOM Polyclonal Antibody	A62158	Human	ELISA, IHC
FUS Polyclonal Antibody	A51449	Human	ELISA, WB, IHC, IF
Fusion Glycoprotein F0 Polyclonal Antibody	A56234	Canine distemper virus	ELISA
Fusion Glycoprotein F0 Polyclonal Antibody	A64438	Human respiratory syncytial virus A	ELISA
Fxn Polyclonal Antibody	A59162	Mouse, Human	ELISA, WB
FXYD1 Polyclonal Antibody	A62618	Human	ELISA
FXYD6 Polyclonal Antibody	A59166	Human	ELISA, IHC
FXYD7 Polyclonal Antibody	A59170	Human	ELISA, IHC, IF
FYCO1 Polyclonal Antibody	A59174	Human, Mouse	ELISA, WB, IHC
FYN Polyclonal Antibody	A59178	Human	ELISA, IHC
FYTTD1 Polyclonal Antibody	A59182	Human	ELISA, WB
FZR1 Polyclonal Antibody	A51405	Human	ELISA, WB, IHC
G0S2 Polyclonal Antibody	A50229	Human, Mouse	ELISA, WB, IHC
G2E3 Polyclonal Antibody	A69931	Human	ELISA, IF
G6PC Polyclonal Antibody	A62622	Human	ELISA, IHC
GABARAP Polyclonal Antibody	A51797	Human	ELISA, WB
GABARAPL2 Polyclonal Antibody	A51745	Human	ELISA, WB
GABARAPL2 Polyclonal Antibody	A51749	Human	ELISA, WB
GABRA4 Polyclonal Antibody	A50233	Human, Mouse	ELISA, WB, IHC
GABRA6 Polyclonal Antibody	A64826	Human	ELISA, IHC
GABRB2 Polyclonal Antibody	A59186	Human	ELISA, IHC, IF
GABRQ Polyclonal Antibody	A68058	Human	ELISA, IHC, IF
Gad2 Polyclonal Antibody	A56935	Mouse	ELISA
gadA Polyclonal Antibody	A53448	Escherichia coli	ELISA
GADD45A Polyclonal Antibody	A56859	Human	ELISA
GADD45B Polyclonal Antibody	A62626	Human	ELISA, IHC
GADD45G Polyclonal Antibody	A57309	Human	ELISA, IHC
gag Polyclonal Antibody	A50708	Human immunodeficiency virus type 1 group M subtype B	ELISA
gag Polyclonal Antibody	A55695	Moloney murine leukemia virus	ELISA

Product Name	Cat. #	Species Reactivity	Applications
gag Polyclonal Antibody	A64434	Moloney murine leukemia virus	ELISA
GAGE5 Polyclonal Antibody	A55403	Human	ELISA, IHC
gag-pol Polyclonal Antibody	A68246	Human immunodeficiency virus type 1 group M subtype B	ELISA
GALE Polyclonal Antibody	A56731	Human	ELISA, WB, IHC, IF
GALK2 Polyclonal Antibody	A59190	Human	ELISA, IHC
GALM Polyclonal Antibody	A55591	Human, Mouse, Rat	ELISA, WB, IHC
GALNS Polyclonal Antibody	A53750	Human	ELISA
GALNT12 Polyclonal Antibody	A59194	Human	ELISA, IHC
GALNT14 Polyclonal Antibody	A51073	Human, Mouse, Rat	ELISA, WB, IHC
GALNT15 Polyclonal Antibody	A66358	Human, Mouse	ELISA, WB, IHC
GALNT16 Polyclonal Antibody	A66946	Human, Mouse	ELISA, WB
GALNT7 Polyclonal Antibody	A65866	Human, Mouse	ELISA, WB, IHC
GALT Polyclonal Antibody	A54772	Human, Mouse, Rat	ELISA, WB, IHC, IF
Gamt Polyclonal Antibody	A54515	Rat, Mouse	ELISA, WB
GAPDH Polyclonal Antibody	A52409	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC, IF
GAPDH Polyclonal Antibody	A52540	Chicken	ELISA
GAR1 Polyclonal Antibody	A67630	Human	ELISA, IHC
GAREM1 Polyclonal Antibody	A67282	Human	ELISA, IHC, IF
GAS2 Polyclonal Antibody	A62630	Human	ELISA
gata1-a Polyclonal Antibody	A68807	Xenopus laevis	ELISA
GATAD2A Polyclonal Antibody	A69019	Human, Rat	ELISA, WB, IHC
GBP4 Polyclonal Antibody	A69131	Human	ELISA, WB, IHC, IF
GBP5 Polyclonal Antibody	A59198	Human, Mouse	ELISA, WB, IHC
GBP6 Polyclonal Antibody	A65534	Human	ELISA, IHC
GBX2 Polyclonal Antibody	A69507	Human	ELISA, IHC
GC Polyclonal Antibody	A57866	Human	ELISA, WB, IHC
GCDH Polyclonal Antibody	A66898	Human, Mouse	ELISA, WB, IHC, IP
GCKR Polyclonal Antibody	A56334	Human	ELISA, IHC
GCNT2 Polyclonal Antibody	A62634	Human	ELISA
GCSAM Polyclonal Antibody	A59202	Human	ELISA, IHC
GCSAML Polyclonal Antibody	A65094	Human	ELISA, IHC
GCY1 Polyclonal Antibody	A50752	Saccharomyces cerevisiae	ELISA
GCY1 Polyclonal Antibody	A50756	Saccharomyces cerevisiae	ELISA
gD Polyclonal Antibody	A55391	Human herpesvirus 1	ELISA
GDAP1L1 Polyclonal Antibody	A59206	Human	ELISA, WB, IHC
GDAP2 Polyclonal Antibody	A51285	Human, Mouse	ELISA, WB
GDF15 Polyclonal Antibody	A52186	Human	ELISA, WB, IHC
GDF2 Polyclonal Antibody	A59210	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
Gdf5 Polyclonal Antibody	A52760	Mouse, Human	ELISA, IHC
GDF9 Polyclonal Antibody	A55170	Human	ELISA, IHC
GDI2 Polyclonal Antibody	A50237	Human, Mouse	ELISA, WB, IHC
GDI2 Polyclonal Antibody	A51496	Human	ELISA, WB
GDPD2 Polyclonal Antibody	A59214	Human	ELISA, WB, IHC, IF
GEMIN4 Polyclonal Antibody	A69511	Human	ELISA, IHC, IF
GEMIN6 Polyclonal Antibody	A66430	Human	ELISA, IHC
GEMIN7 Polyclonal Antibody	A54588	Human	ELISA, WB
GEMIN8 Polyclonal Antibody	A59218	Human, Mouse	ELISA, WB, IHC
Genome Polyprotein Polyclonal Antibody	A56130	Dengue virus	ELISA
Genome Polyprotein Polyclonal Antibody	A56743	Hepatitis C virus genotype 1a	ELISA
Genome Polyprotein Polyclonal Antibody	A68253	Human enterovirus 71	ELISA
Genome Polyprotein Polyclonal Antibody	A68254	Human rhinovirus A serotype 89	ELISA
GFAP Polyclonal Antibody	A52933	Human, Mouse, Rat	ELISA, WB, IHC, IF
GFOD2 Polyclonal Antibody	A50949	Human, Rat	ELISA, WB, IHC
GFRAL Polyclonal Antibody	A65526	Human, Mouse	ELISA, WB, IHC
gG Polyclonal Antibody	A54852	Human herpesvirus 1	ELISA
GGPS1 Polyclonal Antibody	A50241	Human	ELISA, IHC
GGTLC1 Polyclonal Antibody	A62638	Human	ELISA, IHC
GGTLC2 Polyclonal Antibody	A62642	Human, Mouse	ELISA, WB
gH Polyclonal Antibody	A56591	Human cytomegalovirus	ELISA
GH2 Polyclonal Antibody	A59222	Human, Mouse	ELISA, WB, IHC
GHITM Polyclonal Antibody	A59226	Human, Mouse	ELISA, WB, IHC, IF
GHRH Polyclonal Antibody	A50245	Human	ELISA
GHRHR Polyclonal Antibody	A68535	Human	ELISA, IHC, IF
GID8 Polyclonal Antibody	A67746	Human	ELISA, IHC
GIF Polyclonal Antibody	A50249	Human, Mouse	ELISA, WB, IHC
GIF Polyclonal Antibody	A56070	Human	ELISA, IHC
GIMAP7 Polyclonal Antibody	A51049	Human	ELISA, WB
GINS1 Polyclonal Antibody	A69863	Human	ELISA, IHC
GINS2 Polyclonal Antibody	A51409	Human	ELISA, IHC
GINS4 Polyclonal Antibody	A67438	Human	ELISA, IHC
GIT2 Polyclonal Antibody	A64646	Human	ELISA, IHC
GJA4 Polyclonal Antibody	A59230	Human	ELISA, IHC, IF
GJB3 Polyclonal Antibody	A59234	Human, Mouse	ELISA, WB, IHC, IF
GJB4 Polyclonal Antibody	A67306	Human	ELISA
GJB5 Polyclonal Antibody	A50253	Human	ELISA, IHC, IF
GJB6 Polyclonal Antibody	A59238	Human	ELISA, IHC, IF
GJB7 Polyclonal Antibody	A65614	Human	ELISA, IF
GJD2 Polyclonal Antibody	A69307	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
GK2 Polyclonal Antibody	A50257	Human	ELISA, WB, IHC
GK3P Polyclonal Antibody	A50901	Human	ELISA, WB, IHC
GK5 Polyclonal Antibody	A59242	Human	ELISA, WB
GKAP1 Polyclonal Antibody	A59246	Human	ELISA, IHC
GKN1 Polyclonal Antibody	A64128	Human	ELISA
Gkn2 Polyclonal Antibody	A67358	Mouse	ELISA
gL Polyclonal Antibody	A55914	Human cytomegalovirus	ELISA
GLA Polyclonal Antibody	A53220	Human	ELISA, IHC
glaB1 Polyclonal Antibody	A57445	Bacteroides vulgatus	ELISA
GLB1L2 Polyclonal Antibody	A66082	Human	ELISA, IHC
GLCE Polyclonal Antibody	A55279	Human, Mouse, Rat	ELISA, WB, IHC
GLE1 Polyclonal Antibody	A65106	Human	ELISA, WB, IHC, IF
GLI1 Polyclonal Antibody	A52884	Human	ELISA
GLIPR1 Polyclonal Antibody	A52864	Human	ELISA, IF
GLIPR2 Polyclonal Antibody	A67702	Human, Mouse	ELISA, WB
GLP1R Polyclonal Antibody	A53356	Human	ELISA, WB, IHC, IF
GLP1R Polyclonal Antibody	A55798	Human, Mouse	ELISA, WB, IF
glpD Polyclonal Antibody	A53840	Escherichia coli	ELISA
glpK Polyclonal Antibody	A52504	Escherichia coli	ELISA, WB
GLRX Polyclonal Antibody	A57736	Human	ELISA, IHC, IF
GLS Polyclonal Antibody	A62646	Human, Mouse	ELISA, WB, IHC
GLS Polyclonal Antibody	A62650	Human	ELISA
GLS2 Polyclonal Antibody	A67898	Human	ELISA
glsA1 Polyclonal Antibody	A53703	Escherichia coli	ELISA, WB
GLT1D1 Polyclonal Antibody	A66714	Human	ELISA, IHC
GLT8D1 Polyclonal Antibody	A65170	Human, Mouse	ELISA, WB, IHC, IF
GLT8D2 Polyclonal Antibody	A57894	Human, Mouse	ELISA, WB
GLTP Polyclonal Antibody	A62654	Human	ELISA
gluD Polyclonal Antibody	A56843	Clostridium difficile	ELISA
Glycoprotein G Polyclonal Antibody	A53464	Rabies virus	ELISA, WB
Glycoprotein Polyclonal Antibody	A64554	Rabies virus	ELISA
GM2A Polyclonal Antibody	A51505	Human, Rat	ELISA, WB
GMDS Polyclonal Antibody	A69515	Human	ELISA, IHC, IF
GMEB1 Polyclonal Antibody	A59250	Human	ELISA
GMFG Polyclonal Antibody	A62658	Human	ELISA
GMNN Polyclonal Antibody	A59254	Human	ELISA, IHC, IF
GMPPB Polyclonal Antibody	A59258	Human	ELISA, IHC
GMPR Polyclonal Antibody	A55595	Human	ELISA, IHC
GMPR2 Polyclonal Antibody	A51317	Human	ELISA
GNA14 Polyclonal Antibody	A62662	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
GNAI3 Polyclonal Antibody	A62666	Human	ELISA
Gnao1 Polyclonal Antibody	A55105	Mouse, Rat	ELISA, WB
GNAS Polyclonal Antibody	A68803	Human	ELISA, WB, IF
GNAS Polyclonal Antibody	A64422	Human, Mouse, Rat	ELISA, WB, IHC, IF
GNB5 Polyclonal Antibody	A55986	Human	ELISA, IHC, IF
GNG10 Polyclonal Antibody	A59262	Human	ELISA, IHC, IF
GNG11 Polyclonal Antibody	A62670	Human	ELISA, IHC, IF
GNG4 Polyclonal Antibody	A59266	Human	ELISA, IHC, IF
GNG5 Polyclonal Antibody	A62674	Human	ELISA, IHC, IF
GNG7 Polyclonal Antibody	A62678	Human	ELISA
GNGT2 Polyclonal Antibody	A59270	Human	ELISA
GNL1 Polyclonal Antibody	A62682	Human	ELISA, WB, IF
GNL2 Polyclonal Antibody	A57329	Human	ELISA
GNL3L Polyclonal Antibody	A51257	Human	ELISA
GNMT Polyclonal Antibody	A62686	Human	ELISA, IHC, IF
GNPNAT1 Polyclonal Antibody	A67098	Human, Mouse	ELISA, WB, IHC
GNRH2 Polyclonal Antibody	A56278	Human	ELISA, IHC
Goat Serum Albumin Polyclonal Antibody	A57551	Goat	ELISA
GOLM1 Polyclonal Antibody	A54121	Human, Mouse	ELISA, WB, IHC
GOLPH3L Polyclonal Antibody	A59274	Human	ELISA, IHC
GOSR2 Polyclonal Antibody	A50261	Human	ELISA, WB, IHC
GOT1 Polyclonal Antibody	A52800	Human, Mouse	ELISA, WB, IHC, IF
GOT2 Polyclonal Antibody	A51441	Human	ELISA, WB, IHC, IF
GP Polyclonal Antibody	A64246	Zaire ebolavirus	ELISA
GP Polyclonal Antibody	A64254	Zaire ebolavirus	ELISA
GP Polyclonal Antibody	A64414	Zaire ebolavirus	ELISA
Gp2 Polyclonal Antibody	A67678	Mouse	ELISA
GPA33 Polyclonal Antibody	A51213	Human	ELISA, WB, IHC
GPC Polyclonal Antibody	A64474	Lymphocytic choriomeningitis virus	ELISA
GPC4 Polyclonal Antibody	A55239	Mouse	ELISA
GPC4 Polyclonal Antibody	A59278	Human, Mouse	ELISA, WB, IHC, IF
GPC6 Polyclonal Antibody	A55874	Human	ELISA
GPD1 Polyclonal Antibody	A55994	Human	ELISA, IHC, IF
GPD2 Polyclonal Antibody	A62694	Human	ELISA, IHC
GPI Polyclonal Antibody	A53853	Human, Mouse, Rat	ELISA, WB, IHC, IF, IP
Gpi Polyclonal Antibody	A57025	Mouse	ELISA
GPM6A Polyclonal Antibody	A59282	Human, Mouse	ELISA, WB, IF
GPM6B Polyclonal Antibody	A64670	Human, Mouse	ELISA, WB, IHC, IF
GPR1 Polyclonal Antibody	A68543	Human	ELISA, IHC, IF
GPR101 Polyclonal Antibody	A66822	Human	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
GPR141 Polyclonal Antibody	A65906	Human	ELISA
GPR146 Polyclonal Antibody	A59286	Human	ELISA, IHC, IF
GPR156 Polyclonal Antibody	A66778	Human	ELISA, IHC
GPR161 Polyclonal Antibody	A53973	Human	ELISA, WB, IF
GPR17 Polyclonal Antibody	A59290	Human, Mouse	ELISA, WB, IF
GPR173 Polyclonal Antibody	A67734	Human	ELISA, IHC, IF
GPR176 Polyclonal Antibody	A64742	Human	ELISA, WB, IF
GPR18 Polyclonal Antibody	A59294	Human, Mouse	ELISA, WB, IHC, IF
GPR183 Polyclonal Antibody	A62698	Human	ELISA, IHC, IF
GPR21 Polyclonal Antibody	A67162	Human	ELISA, IHC, IF
GPR3 Polyclonal Antibody	A57914	Human	ELISA, IHC, IF
GPR34 Polyclonal Antibody	A68062	Human, Mouse	ELISA, WB, IF
GPR75 Polyclonal Antibody	A59298	Human	ELISA, WB, IHC, IF
GPR82 Polyclonal Antibody	A67122	Human, Mouse	ELISA, WB, IHC, IF
GPR83 Polyclonal Antibody	A67330	Human	ELISA, WB, IF
GPR84 Polyclonal Antibody	A67298	Human	ELISA, IF
GPRC5B Polyclonal Antibody	A59302	Human	ELISA, IF
GPSM1 Polyclonal Antibody	A65966	Human	ELISA
GPSM3 Polyclonal Antibody	A68158	Human	ELISA, IHC
GPT Polyclonal Antibody	A56418	Human, Rat	ELISA, WB, IHC, IF
GPX7 Polyclonal Antibody	A59306	Human	ELISA, IHC
GPX8 Polyclonal Antibody	A66886	Human	ELISA, IHC, IF
GRA1 Polyclonal Antibody	A55806	Toxoplasma gondii	ELISA
GRAMD1B Polyclonal Antibody	A64934	Human	ELISA, IHC, IF
GRB2 Polyclonal Antibody	A62702	Human	ELISA
GREM1 Polyclonal Antibody	A69519	Human	ELISA, IHC
GRHL2 Polyclonal Antibody	A65506	Human	ELISA, IHC, IF
GRIA3 Polyclonal Antibody	A50265	Human	ELISA, IHC, IF
GRIN2A Polyclonal Antibody	A52916	Human, Mouse, Rat	ELISA, WB, IHC, IF
GRK4 Polyclonal Antibody	A50269	Human	ELISA, WB, IHC, IF
groL Polyclonal Antibody	A50857	Yersinia pestis	ELISA
GRP Polyclonal Antibody	A54404	Human	ELISA, IHC, IF
GRP Polyclonal Antibody	A57181	Human	ELISA, IHC
GRPR Polyclonal Antibody	A68547	Human	ELISA, IF
GSC Polyclonal Antibody	A69523	Human	ELISA, IHC
GSDMB Polyclonal Antibody	A59314	Human	ELISA, WB, IHC
GSDMC Polyclonal Antibody	A67802	Human	ELISA, IHC
GSDMD Polyclonal Antibody	A59318	Human	ELISA, WB
gshA Polyclonal Antibody	A64522	Escherichia coli	ELISA
GSK3B Polyclonal Antibody	A62706	Human, Mouse	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
GSN Polyclonal Antibody	A50273	Human	ELISA, WB, IHC, IF
GSS Polyclonal Antibody	A69527	Human, Mouse, Rat	ELISA, WB, IHC
GSTA1 Polyclonal Antibody	A62710	Human	ELISA, IHC, IF
GSTA4 Polyclonal Antibody	A50277	Human	ELISA, IHC, IF
GSTK1 Polyclonal Antibody	A57249	Human	ELISA, WB
Gstp1 Polyclonal Antibody	A52784	Rat, Human	ELISA, WB
GSTP1 Polyclonal Antibody	A54109	Human	ELISA, WB
GSTP1 Polyclonal Antibody	A54113	Human	ELISA
Gstp1 Polyclonal Antibody	A56198	Mouse, Human	ELISA, WB
GSTT1 Polyclonal Antibody	A59322	Human	ELISA, IHC, IF
GT-1 Polyclonal Antibody	A54882	Arabidopsis thaliana	ELISA
GTDC1 Polyclonal Antibody	A65018	Human	ELISA, WB
GTF2B Polyclonal Antibody	A54471	Human	ELISA, WB, IHC, IF
GTF2E2 Polyclonal Antibody	A53643	Human, Mouse	ELISA, WB
GTF2F2 Polyclonal Antibody	A62714	Human	ELISA, WB
GTF2H2 Polyclonal Antibody	A64734	Human	ELISA
GTF2H5 Polyclonal Antibody	A54684	Human	ELISA, IHC, IF
GTF3C5 Polyclonal Antibody	A69331	Human	ELISA, IHC
GTF3C6 Polyclonal Antibody	A59326	Human	ELISA, IHC
GTPBP1 Polyclonal Antibody	A69531	Human	ELISA, IHC, IF
GTPBP6 Polyclonal Antibody	A59330	Human	ELISA, WB, IHC
GTSF1 Polyclonal Antibody	A66890	Human	ELISA
GUCA1A Polyclonal Antibody	A62718	Human	ELISA
Guca2b Polyclonal Antibody	A56330	Mouse	ELISA
GUCA2B Polyclonal Antibody	A56342	Human	ELISA, IHC
Guinea Pig Hemoglobin Polyclonal Antibody	A57569	Guinea Pig	ELISA, WB
Guinea Pig Serum Albumin Polyclonal Antibody	A57556	Guinea Pig	ELISA
GV51_0914 Polyclonal Antibody	A64334	Gardnerella vaginalis 5-1	ELISA
GYPA Polyclonal Antibody	A53352	Human	ELISA
GYPE Polyclonal Antibody	A62722	Human	ELISA
Gzma Polyclonal Antibody	A59334	Mouse	ELISA
GZMB Polyclonal Antibody	A55359	Human, Mouse	ELISA, WB
H2AFB1 Polyclonal Antibody	A62726	Human	ELISA, IHC, IF
H2AFV Polyclonal Antibody	A64060	Human	ELISA
H2AFY2 Polyclonal Antibody	A67878	Human	ELISA
H2-D1 Polyclonal Antibody	A57413	Mouse, Human	ELISA, WB
H2-K1 Polyclonal Antibody	A53742	Mouse	ELISA
HA Polyclonal Antibody	A55515	Influenza A virus (H7N9)	ELISA
HA Polyclonal Antibody	A64446	Influenza A virus	ELISA
HABP2 Polyclonal Antibody	A50688	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
HACD2 Polyclonal Antibody	A65662	Human	ELISA, WB, IHC, IF
HACE1 Polyclonal Antibody	A69055	Human	ELISA, IHC
HADHB Polyclonal Antibody	A51593	Human	ELISA, IHC
HAGH Polyclonal Antibody	A51970	Human	ELISA, IHC
HAL Polyclonal Antibody	A69535	Human	ELISA, IHC, IF
HAO1 Polyclonal Antibody	A69311	Human	ELISA, IHC, IF
HARS2 Polyclonal Antibody	A54592	Human, Mouse	ELISA, WB, IHC
HAS1 Polyclonal Antibody	A68551	Human	ELISA, IF
Has2 Polyclonal Antibody	A55086	Mouse	ELISA
HAVCR1 Polyclonal Antibody	A50652	Human, Mouse	ELISA, WB, IHC
Havcr1 Polyclonal Antibody	A52368	Mouse, Human	ELISA, IHC
HAVCR2 Polyclonal Antibody	A52552	Human, Mouse	ELISA, IHC, IF
HAX1 Polyclonal Antibody	A51721	Human	ELISA, WB, IHC, IF
Hba Polyclonal Antibody	A53679	Mouse, Human	ELISA, WB
HBA1 Polyclonal Antibody	A57842	Human	ELISA
HBB Polyclonal Antibody	A50281	Human	ELISA, IHC
Hbb-b2 Polyclonal Antibody	A53598	Mouse	ELISA
HBD Polyclonal Antibody	A59338	Human	ELISA, IHC
HBE1 Polyclonal Antibody	A59342	Human	ELISA, IHC
HBEGF Polyclonal Antibody	A69147	Human	ELISA, IF
HBG1 Polyclonal Antibody	A55715	Human	ELISA, IHC
HBQ1 Polyclonal Antibody	A59346	Human	ELISA
HCFC1R1 Polyclonal Antibody	A67322	Human	ELISA, IHC
HDAC1 Polyclonal Antibody	A68304	Human, Mouse, Zebrafish	ELISA, WB, IHC, ChIP
HDAC3 Polyclonal Antibody	A68305	Human, Mouse	ELISA, WB, IHC, IF, ChIP
HDAC4 Polyclonal Antibody	A52524	Human	ELISA, IHC
HDAC7 Polyclonal Antibody	A52356	Human	ELISA
HDAC8 Polyclonal Antibody	A55962	Human	ELISA, IHC, IF
HDDC3 Polyclonal Antibody	A51085	Human, Mouse	ELISA, WB, IHC
HDGF Polyclonal Antibody	A53048	Human	ELISA
HDHD3 Polyclonal Antibody	A59354	Human, Mouse	ELISA, WB, IHC
HEATR9 Polyclonal Antibody	A59358	Human	ELISA, WB, IHC
HEBP2 Polyclonal Antibody	A57177	Human	ELISA, IHC
HECTD2 Polyclonal Antibody	A69923	Human	ELISA, IHC, IF
HEMK1 Polyclonal Antibody	A59362	Human	ELISA, IHC
HEPACAM Polyclonal Antibody	A68555	Human	ELISA, WB, IF
HEPACAM2 Polyclonal Antibody	A62730	Human	ELISA, IHC
HERC3 Polyclonal Antibody	A64746	Human	ELISA
HERC6 Polyclonal Antibody	A66094	Human	ELISA, IHC
HESX1 Polyclonal Antibody	A67430	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
HEXB Polyclonal Antibody	A68559	Human	ELISA, IHC
HEXIM2 Polyclonal Antibody	A62734	Human, Mouse, Rat	ELISA, WB, IHC
HEY1 Polyclonal Antibody	A69347	Human	ELISA, IHC
HFE Polyclonal Antibody	A64894	Human	ELISA
HHIPL1 Polyclonal Antibody	A66306	Human	ELISA
HHLA3 Polyclonal Antibody	A59366	Human	ELISA
HIF1A Polyclonal Antibody	A56651	Human	ELISA, WB, IHC, IF
HIF1AN Polyclonal Antibody	A59374	Human, Mouse	ELISA, WB, IHC, IF, IP
HIGD1B Polyclonal Antibody	A68006	Human	ELISA, IHC, IF
HINFP Polyclonal Antibody	A69539	Human, Rat	ELISA, WB
HIP1R Polyclonal Antibody	A68563	Human, Rat	ELISA, WB, IHC
HIPK1 Polyclonal Antibody	A69027	Human	ELISA, IHC, IF
HIST1H2AG Polyclonal Antibody	A68306	Human	ELISA, WB, IHC, IF, ChIP
HIST1H2BK Polyclonal Antibody	A62738	Human	ELISA, IHC
HIST2H2AC Polyclonal Antibody	A62742	Human	ELISA, IHC
Histidine-Rich Protein PFHRP-II Polyclonal Antibody	A64450	Plasmodium falciparum	ELISA
Histone Acetyltransferase KAT5 Polyclonal Antibody	A54421	Human	ELISA, IHC, IF
Histone Acetyltransferase KAT7 Polyclonal Antibody	A56639	Human, Rat	ELISA, WB, IHC
Histone H2A.J Polyclonal Antibody	A53292	Human, Mouse	ELISA, WB, IHC
Histone H2A.x Polyclonal Antibody	A54189	Human	ELISA, IHC
Histone H2B Polyclonal Antibody	A57193	Cairina moschata	ELISA
Histone H2B Type 1-B Polyclonal Antibody	A50289	Human	ELISA, IHC, IF
Histone H2B Type 1-C/E/F/G/I Polyclonal Antibody	A51588	Human	ELISA, WB
Histone H2B Type 1-M Polyclonal Antibody	A52528	Mouse	ELISA
Histone H2B Type 1-N Polyclonal Antibody	A51217	Human	ELISA, IHC
Histone H3.3 Polyclonal Antibody	A52644	Human	ELISA, WB, IHC
Histone H3.3 Polyclonal Antibody	A55567	Human	ELISA, WB, IHC
HJURP Polyclonal Antibody	A56118	Human	ELISA, WB
HK1 Polyclonal Antibody	A54337	Human	ELISA
HK1 Polyclonal Antibody	A54597	Human, Rat	ELISA, WB
HKDC1 Polyclonal Antibody	A64866	Human	ELISA, WB, IHC
HLA-B Polyclonal Antibody	A52237	Human	ELISA, WB, IHC, IF
HLA-C Polyclonal Antibody	A52484	Human	ELISA, WB, IHC, IF
HLA-DMA Polyclonal Antibody	A59378	Human	ELISA, IHC
HLA-DPA1 Polyclonal Antibody	A53174	Human	ELISA, WB, IF
HLA-DPB1 Polyclonal Antibody	A57886	Human	ELISA, WB, IHC
HLA-DQA1 Polyclonal Antibody	A52488	Human	ELISA, IHC
HLA-DQA2 Polyclonal Antibody	A52480	Human	ELISA, IHC
HLA-DRA Polyclonal Antibody	A53149	Human	ELISA, IHC
HLA-DRB1 Polyclonal Antibody	A50716	Human	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
HLA-DRB1 Polyclonal Antibody	A52026	Human	ELISA
HLA-DRB4 Polyclonal Antibody	A59382	Human	ELISA, WB, IHC
HLA-E Polyclonal Antibody	A50748	Human	ELISA
HLA-G Polyclonal Antibody	A53328	Human	ELISA, WB, IHC, IF
hly Polyclonal Antibody	A56242	Staphylococcus aureus	ELISA
HM13 Polyclonal Antibody	A59386	Human, Mouse	ELISA, WB, IF
HMGB1 Polyclonal Antibody	A51497	Human, Mouse	ELISA, WB, IHC
HMGB4 Polyclonal Antibody	A51173	Human	ELISA, IHC
HMGCLL1 Polyclonal Antibody	A59390	Human	ELISA
HMGCR Polyclonal Antibody	A52932	Human	ELISA, WB, IHC, IF
Hmgcs1 Polyclonal Antibody	A51029	Mouse	ELISA
Hmgcs1 Polyclonal Antibody	A51033	Mouse	ELISA, WB
HMGCS2 Polyclonal Antibody	A68567	Human, Rat	ELISA, WB, IHC, IF
HMGN3 Polyclonal Antibody	A59394	Human	ELISA, IHC
HMOX1 Polyclonal Antibody	A52544	Human	ELISA, IHC, IF
HN1 Polyclonal Antibody	A56535	Human, Rat	ELISA, WB
HNRNPA1 Polyclonal Antibody	A53428	Human, Mouse	ELISA, WB, IHC, IF
HNRNPA2B1 Polyclonal Antibody	A54605	Human	ELISA, IHC, IF
HNRNPAB Polyclonal Antibody	A69543	Human	ELISA, WB, IHC, IF
HNRNPD Polyclonal Antibody	A53913	Human	ELISA, WB, IHC
HNRNPH1 Polyclonal Antibody	A51825	Human	ELISA, WB
HNRNPH1 Polyclonal Antibody	A56615	Human	ELISA, WB, IHC
HNRNPL Polyclonal Antibody	A50293	Human	ELISA, IHC, IF
HNRNPL Polyclonal Antibody	A59398	Human, Mouse	ELISA, WB, IHC
HOMER1 Polyclonal Antibody	A52492	Human	ELISA, IHC
HOMER2 Polyclonal Antibody	A59402	Human	ELISA, IHC
HOMER3 Polyclonal Antibody	A59406	Human, Mouse	ELISA, WB, IHC
HOOK1 Polyclonal Antibody	A68046	Human	ELISA, WB, IHC
HOPX Polyclonal Antibody	A59410	Human	ELISA, IHC
HORMAD1 Polyclonal Antibody	A65998	Human	ELISA
HORMAD2 Polyclonal Antibody	A66494	Human	ELISA, IHC
Horse Serum Albumin Polyclonal Antibody	A57536	Horse	ELISA
HOXA4 Polyclonal Antibody	A68571	Human, Rat	ELISA, WB, IHC
HOXA6 Polyclonal Antibody	A50297	Human, Mouse	ELISA, WB, IHC
HOXB4 Polyclonal Antibody	A68575	Human	ELISA, IHC, IF
HOXC11 Polyclonal Antibody	A59414	Human	ELISA, IHC, IF
HOXC9 Polyclonal Antibody	A62746	Human	ELISA
HOXD1 Polyclonal Antibody	A70311	Human, Mouse, Rat	ELISA, WB, IHC
HOXD3 Polyclonal Antibody	A68579	Human, Rat	ELISA, WB
HOXD8 Polyclonal Antibody	A68583	Human, Mouse	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
Hp Polyclonal Antibody	A52142	Rat, Mouse	ELISA, WB, IHC
HP Polyclonal Antibody	A55503	Human, Mouse	ELISA, WB, IHC
HPCAL1 Polyclonal Antibody	A59418	Human	ELISA, IHC, IF
HPCAL4 Polyclonal Antibody	A68082	Human, Mouse	ELISA, WB
HPD Polyclonal Antibody	A52564	Human, Mouse	ELISA, WB, IHC, IF
HPN Polyclonal Antibody	A69547	Human, Mouse, Rat	ELISA, WB, IHC, IF
HPRT1 Polyclonal Antibody	A52336	Human	ELISA, IHC
HPS6 Polyclonal Antibody	A69023	Human	ELISA, IHC, IF
Hpx Polyclonal Antibody	A52880	Human, Rat	ELISA, WB
Hras Polyclonal Antibody	A54527	Rat, Mouse, Human	ELISA, WB, IHC
HRSP12 Polyclonal Antibody	A62750	Human	ELISA, IHC, IF
HS2ST1 Polyclonal Antibody	A59422	Human, Mouse, Rat	ELISA, WB, IHC
HS3ST1 Polyclonal Antibody	A62754	Human	ELISA, IHC
HSBP1 Polyclonal Antibody	A53001	Human	ELISA, IHC
HSD11B1 Polyclonal Antibody	A56290	Human	ELISA, WB, IHC, IF
HSD11B1L Polyclonal Antibody	A65834	Human	ELISA, IHC
HSD17B10 Polyclonal Antibody	A56106	Human, Mouse	ELISA, WB
HSD17B11 Polyclonal Antibody	A56751	Human, Mouse	ELISA, WB
HSD17B12 Polyclonal Antibody	A59426	Human	ELISA, WB, IHC, IF
HSD17B14 Polyclonal Antibody	A57273	Human, Mouse	ELISA, WB
HSDL2 Polyclonal Antibody	A59430	Human	ELISA, WB, IHC
HSP90AB1 Polyclonal Antibody	A50001	Human, Mouse, Rat, Zebrafish	ELISA, WB
HSP90B1 Polyclonal Antibody	A53162	Human	ELISA, WB, IHC
HSPA12B Polyclonal Antibody	A64132	Human	ELISA, IHC
HSPA13 Polyclonal Antibody	A54141	Human	ELISA, IHC
Hspa1b Polyclonal Antibody	A59434	Mouse, Human	ELISA, WB
HSPA5 Polyclonal Antibody	A53981	Human	ELISA, IHC
HSPA6 Polyclonal Antibody	A53804	Human	ELISA, WB
HSPB7 Polyclonal Antibody	A59438	Human	ELISA, IHC
HSPBAP1 Polyclonal Antibody	A51177	Human	ELISA, IHC
HSPE1 Polyclonal Antibody	A51986	Human	ELISA
HTR1E Polyclonal Antibody	A62758	Human	ELISA, WB, IHC, IF
HTR2A Polyclonal Antibody	A50301	Human	ELISA, IHC, IF
HTR3D Polyclonal Antibody	A65490	Human	ELISA, WB, IF
Human Serum Albumin Polyclonal Antibody	A57544	Human	ELISA
Human Serum Albumin Polyclonal Antibody	A57748	Human	ELISA
hup1 Polyclonal Antibody	A50829	Streptomyces lividans	ELISA
HUS1 Polyclonal Antibody	A54483	Human	ELISA, WB, IHC, IF
HYDIN Polyclonal Antibody	A59442	Human	ELISA
HYPM Polyclonal Antibody	A62358	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
IAPP Polyclonal Antibody	A52241	Human, Mouse	ELISA, WB, IHC
ICA1 Polyclonal Antibody	A53268	Human, Mouse	ELISA, WB, IHC
ICA1L Polyclonal Antibody	A59446	Human	ELISA, WB
ICT1 Polyclonal Antibody	A55587	Human	ELISA, WB, IHC
ID2 Polyclonal Antibody	A54860	Human	ELISA, IHC, IF
IDH1 Polyclonal Antibody	A53345	Mouse, Human	ELISA, IHC, IF
IDH3B Polyclonal Antibody	A53686	Human, Mouse	ELISA, WB
IDH3G Polyclonal Antibody	A56026	Human	ELISA
IDI2 Polyclonal Antibody	A59450	Human	ELISA, IHC
IDO1 Polyclonal Antibody	A53757	Human	ELISA, WB, IHC, IF
IDO2 Polyclonal Antibody	A56410	Human	ELISA, IHC, IF
IDUA Polyclonal Antibody	A56066	Human	ELISA, IHC
IER2 Polyclonal Antibody	A59454	Human	ELISA
IFI27L2 Polyclonal Antibody	A59458	Human	ELISA, IHC, IF
IFI44 Polyclonal Antibody	A62762	Human, Mouse	ELISA, WB, IHC, IF
IFIH1 Polyclonal Antibody	A67670	Human	ELISA, IHC, IF
IFIT3 Polyclonal Antibody	A55830	Human	ELISA, WB, IHC, IF
IFIT5 Polyclonal Antibody	A59462	Human	ELISA, WB, IHC, IF
IFITM1 Polyclonal Antibody	A51681	Human	ELISA, WB, IF
IFITM2 Polyclonal Antibody	A62766	Human	ELISA, WB, IHC, IF
IFNA1 Polyclonal Antibody	A56811	Human	ELISA
IFNA10 Polyclonal Antibody	A54990	Human, Mouse	ELISA, WB, IHC
IFNA14 Polyclonal Antibody	A54708	Human, Mouse	ELISA, WB, IHC
IFNA14 Polyclonal Antibody	A62770	Human	ELISA
IFNA2 Polyclonal Antibody	A53753	Human, Mouse	ELISA, WB, IHC, IF
IFNA21 Polyclonal Antibody	A53816	Human	ELISA, WB
IFNA4 Polyclonal Antibody	A56959	Human, Mouse	ELISA, WB, IHC
IFNAG Polyclonal Antibody	A64390	Bovine	ELISA, WB
IFNAR2 Polyclonal Antibody	A50680	Human	ELISA, IHC
IFNB1 Polyclonal Antibody	A51929	Human	ELISA, IHC
Ifne Polyclonal Antibody	A55174	Mouse	ELISA
IFNG Polyclonal Antibody	A51717	Human	ELISA, WB
IFNG Polyclonal Antibody	A51990	Human	ELISA
Ifng Polyclonal Antibody	A52046	Rat	ELISA
IFNG Polyclonal Antibody	A52054	Chicken	ELISA
Ifng Polyclonal Antibody	A52440	Mouse, Human	ELISA, IHC
IFNG Polyclonal Antibody	A53166	Human	ELISA, WB
Ifng Polyclonal Antibody	A57626	Guinea Pig	ELISA, WB
IFNT1 Polyclonal Antibody	A57397	Bovine	ELISA
IFRD2 Polyclonal Antibody	A59466	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
IFT22 Polyclonal Antibody	A51249	Human	ELISA, IHC
IFT27 Polyclonal Antibody	A51321	Human	ELISA, IHC
IFT80 Polyclonal Antibody	A67882	Human	ELISA, IHC
IGF1 Polyclonal Antibody	A52520	Human	ELISA
IGF1R Polyclonal Antibody	A64112	Human	ELISA
Igfbp1 Polyclonal Antibody	A50302	Mouse, Human	ELISA, WB
IGFBP1 Polyclonal Antibody	A54583	Human	ELISA, WB, IHC
igg Heavy Chain Polyclonal Antibody	A57584	goat	ELISA, WB
igg Heavy Chain Polyclonal Antibody	A57789	Human	ELISA
igg Light Chain Polyclonal Antibody	A57596	goat	ELISA
igg Light Chain Polyclonal Antibody	A57793	Human	ELISA, WB
IGHG1 Polyclonal Antibody	A62774	Human	ELISA, IHC
IGLL5 Polyclonal Antibody	A64574	Human	ELISA
IGSF11 Polyclonal Antibody	A59470	Human, Mouse	ELISA, WB
IGSF6 Polyclonal Antibody	A62778	Human	ELISA
IGSF9 Polyclonal Antibody	A69807	Human	ELISA, IHC
IHH Polyclonal Antibody	A54728	Human	ELISA, IHC, IF
III Polyclonal Antibody	A64208	Enterobacteria phage M13	ELISA
IKZF1 Polyclonal Antibody	A64786	Human	ELISA
IL10 Polyclonal Antibody	A56298	Bovine	ELISA
IL10 Polyclonal Antibody	A56495	Macaca mulatta	ELISA
IL12B Polyclonal Antibody	A53304	Sheep	ELISA
IL12B Polyclonal Antibody	A53559	Sheep	ELISA, WB, IHC
IL12B Polyclonal Antibody	A54021	Human, Mouse	ELISA, WB, IHC
IL12RB2 Polyclonal Antibody	A69159	Human	ELISA, IF
IL15 Polyclonal Antibody	A52376	Human	ELISA, IHC
IL18 Polyclonal Antibody	A53635	Human	ELISA, IHC
IL18 Polyclonal Antibody	A54085	Human	ELISA
IL18 Polyclonal Antibody	A56563	Dog	ELISA
IL18BP Polyclonal Antibody	A52126	Human	ELISA, IHC
IL18BP Polyclonal Antibody	A52146	Mouse, Human	ELISA, IHC
IL1A Polyclonal Antibody	A50306	Macaca mulatta	ELISA, WB
IL1A Polyclonal Antibody	A51641	Human	ELISA
IL1B Polyclonal Antibody	A51537	Human	ELISA
Il1b Polyclonal Antibody	A52348	Mouse	ELISA
IL1B Polyclonal Antibody	A53690	Sheep	ELISA, WB
IL1B Polyclonal Antibody	A54680	Sheep	ELISA, WB
IL1F10 Polyclonal Antibody	A56775	Human	ELISA, IHC
IL1RAPL1 Polyclonal Antibody	A69551	Human	ELISA, WB, IHC, IF
IL1RL1 Polyclonal Antibody	A55146	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
IL1RL2 Polyclonal Antibody	A67966	Human	ELISA, IHC
IL1RN Polyclonal Antibody	A52233	Human	ELISA, WB
IL2 Polyclonal Antibody	A59474	Human	ELISA, IHC
IL22RA1 Polyclonal Antibody	A51069	Human	ELISA, IHC
IL24 Polyclonal Antibody	A64634	Human	ELISA
IL2rb Polyclonal Antibody	A54212	Mouse	ELISA
IL2RG Polyclonal Antibody	A50310	Human	ELISA, WB, IHC
IL3 Polyclonal Antibody	A55142	Human	ELISA
IL-33 Polyclonal Antibody	A57721	Sus scrofa, Mouse	ELISA, WB
IL36RN Polyclonal Antibody	A56547	Human, Mouse, Rat	ELISA, WB
IL4 Polyclonal Antibody	A51845	Human	ELISA
IL4 Polyclonal Antibody	A53452	Dog	ELISA
IL6 Polyclonal Antibody	A54025	Human, Mouse	ELISA, WB, IHC
ILF2 Polyclonal Antibody	A52712	Human	ELISA, WB, IHC
ILK Polyclonal Antibody	A52452	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC, IF
ILKAP Polyclonal Antibody	A54596	Human, Mouse, Rat	ELISA, WB
ILVBL Polyclonal Antibody	A62786	Human	ELISA, WB, IHC, IF
Immunoglobulin G2a Polyclonal Antibody	A50221	Mouse	ELISA
IMPACT Polyclonal Antibody	A68014	Human, Mouse	ELISA, WB, IHC
IMPDH2 Polyclonal Antibody	A57797	Human	ELISA, WB, IHC, IF
IMPDH2 Polyclonal Antibody	A59478	Human, Mouse	ELISA, WB, IHC, IF
IMUP Polyclonal Antibody	A53667	Human	ELISA
INCA1 Polyclonal Antibody	A64606	Human	ELISA, IHC
INHA Polyclonal Antibody	A53008	Human	ELISA, IHC
inhA Polyclonal Antibody	A56218	Mycobacterium tuberculosis	ELISA
Inhibin Beta B Chain Polyclonal Antibody	A57644	Human	ELISA, WB, IHC, IF
inlA Polyclonal Antibody	A65630	Listeria monocytogenes serotype 4b	ELISA, WB
INMT Polyclonal Antibody	A53776	Human, Mouse	ELISA, WB, IHC
INPP1 Polyclonal Antibody	A59482	Human	ELISA, IHC
INPP5A Polyclonal Antibody	A55251	Human	ELISA, IHC, IF
INPP5B Polyclonal Antibody	A69555	Human	ELISA, IHC, IF
INS Polyclonal Antibody	A55687	Human	ELISA, IHC
INSL4 Polyclonal Antibody	A64798	Human	ELISA, IHC
Intermediate Capsid Protein VP6 Polyclonal Antibody	A55018	Rotavirus A	ELISA
Intermediate Capsid Protein VP6 Polyclonal Antibody	A56883	Rotavirus A	ELISA
Intermediate Capsid Protein VP6 Polyclonal Antibody	A57101	Rotavirus A	ELISA
INTS6 Polyclonal Antibody	A68587	Human	ELISA, IF
iolG Polyclonal Antibody	A54624	Bacillus subtilis	ELISA, WB
ipaH9.8 Polyclonal Antibody	A55475	Shigella flexneri serotype X	ELISA

Product Name	Cat. #	Species Reactivity	Applications
IPO11 Polyclonal Antibody	A59486	Human	ELISA
IPO13 Polyclonal Antibody	A62790	Human	ELISA, WB, IHC, IF
IQCC Polyclonal Antibody	A65030	Human	ELISA, IHC
IQCF2 Polyclonal Antibody	A66142	Human	ELISA
Irf5 Polyclonal Antibody	A62794	Mouse	ELISA
Irf8 Polyclonal Antibody	A50314	Mouse	ELISA, WB
ISCA2 Polyclonal Antibody	A65798	Human	ELISA, WB, IHC
ISCU Polyclonal Antibody	A67942	Human	ELISA
ISG15 Polyclonal Antibody	A53056	Human	ELISA, WB
IST1 Polyclonal Antibody	A69563	Human	ELISA, IHC, IF
ISX Polyclonal Antibody	A50318	Human	ELISA, IHC
ITGA2 Polyclonal Antibody	A59490	Human	ELISA, IHC
ITGA2B Polyclonal Antibody	A57409	Human	ELISA, IHC
Itgal Polyclonal Antibody	A52868	Mouse	ELISA
ITGAL Polyclonal Antibody	A62798	Human	ELISA, WB, IHC, IF
Itgb1 Polyclonal Antibody	A55150	Rat	ELISA
ITGB1 Polyclonal Antibody	A62802	Mouse, Human	ELISA, WB, IF
ITGB1BP2 Polyclonal Antibody	A68110	Human	ELISA, IHC
ITGB6 Polyclonal Antibody	A68591	Human, Rat	ELISA, WB, IHC, IF
ITIH4 Polyclonal Antibody	A54870	Pig	ELISA
ITIH4 Polyclonal Antibody	A56354	Human	ELISA
ITIH4 Polyclonal Antibody	A56426	Human	ELISA, IHC
ITIH5 Polyclonal Antibody	A59494	Human	ELISA, IHC
ITM2A Polyclonal Antibody	A62806	Human	ELISA, WB, IHC, IF
IZUMO1 Polyclonal Antibody	A66106	Human	ELISA, IHC, IF
IZUMO4 Polyclonal Antibody	A56142	Human, Rat	ELISA, WB, IHC, IF
JAGN1 Polyclonal Antibody	A66190	Human	ELISA, WB, IHC, IF
JAK2 Polyclonal Antibody	A53469	Human	ELISA
JAK3 Polyclonal Antibody	A59498	Human	ELISA
JAM2 Polyclonal Antibody	A52912	Human	ELISA, WB
JPH1 Polyclonal Antibody	A69227	Human	ELISA, IHC, IF
JUP Polyclonal Antibody	A50692	Human	ELISA, WB, IHC
Kallikrein-10 Polyclonal Antibody	A57648	Human	ELISA
KARS Polyclonal Antibody	A64076	Human	ELISA, WB, IHC, IF
KATNAL2 Polyclonal Antibody	A59502	Human	ELISA, WB, IHC
KBTBD2 Polyclonal Antibody	A51025	Human	ELISA, WB
KBTBD4 Polyclonal Antibody	A59506	Human, Mouse	ELISA, WB, IHC
KCMF1 Polyclonal Antibody	A70259	Human	ELISA, IHC
KCNA1 Polyclonal Antibody	A52945	Human	ELISA
KCNA6 Polyclonal Antibody	A62810	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
KCNC3 Polyclonal Antibody	A68875	Human	ELISA, IHC, IF
KCND1 Polyclonal Antibody	A52573	Human	ELISA, IHC, IF
KCNE2 Polyclonal Antibody	A54400	Rat	ELISA
KCNE3 Polyclonal Antibody	A68775	Human	ELISA, IHC
KCNH7 Polyclonal Antibody	A59510	Human	ELISA, IHC, IF
KCNIP4 Polyclonal Antibody	A65522	Human	ELISA, IHC, IF
KCNJ1 Polyclonal Antibody	A53694	Human	ELISA, WB, IF
KCNJ10 Polyclonal Antibody	A56162	Human	ELISA, IHC, IF
KCNJ11 Polyclonal Antibody	A64818	Human, Mouse	ELISA, WB, IHC
KCNJ5 Polyclonal Antibody	A69559	Human, Mouse	ELISA, WB, IHC, IF
KCNJ8 Polyclonal Antibody	A68827	Human	ELISA, IHC, IF
KCNK13 Polyclonal Antibody	A67206	Human	ELISA, WB, IHC, IF
KCNK3 Polyclonal Antibody	A68595	Human	ELISA, IHC
KCNMA1 Polyclonal Antibody	A64666	Human	ELISA, IHC, IF
KCNMB1 Polyclonal Antibody	A59514	Human	ELISA, IF
KCNRG Polyclonal Antibody	A59518	Human	ELISA, IHC
KCNS2 Polyclonal Antibody	A59522	Human, Mouse	ELISA, WB, IF
KCNT1 Polyclonal Antibody	A68911	Human	ELISA, IHC, IF
KCNV2 Polyclonal Antibody	A70119	Human	ELISA, IHC, IF
KCTD1 Polyclonal Antibody	A65446	Human	ELISA
KCTD16 Polyclonal Antibody	A65262	Human	ELISA, IHC
KCTD18 Polyclonal Antibody	A65618	Human	ELISA
KCTD19 Polyclonal Antibody	A64850	Human	ELISA, IHC
KCTD4 Polyclonal Antibody	A59526	Human	ELISA, IHC
KCTD6 Polyclonal Antibody	A66394	Human	ELISA, IHC
KCTD9 Polyclonal Antibody	A65414	Human	ELISA, WB
KDM5A Polyclonal Antibody	A55247	Human	ELISA, IHC, IF
KDM8 Polyclonal Antibody	A59530	Human	ELISA, IHC
KDR Polyclonal Antibody	A55495	Human	ELISA, IHC
KEAP1 Polyclonal Antibody	A50322	Human, Mouse	ELISA, WB, IHC
Keap1 Polyclonal Antibody	A59534	Mouse, Human	ELISA, WB
KEL Polyclonal Antibody	A59538	Human	ELISA, WB, IHC
KHDC1 Polyclonal Antibody	A65138	Human, Mouse	ELISA, WB, IHC, IF
KHDRBS2 Polyclonal Antibody	A65290	Human, Mouse	ELISA, WB, IHC
KIAA0101 Polyclonal Antibody	A54240	Human	ELISA
KIAA0319 Polyclonal Antibody	A68779	Human	ELISA, IHC, IF
KIAA0319L Polyclonal Antibody	A59542	Human	ELISA, IHC, IF
KIAA0513 Polyclonal Antibody	A62814	Human, Mouse	ELISA, WB, IF
KIAA0907 Polyclonal Antibody	A59546	Human, Mouse	ELISA, WB, IHC
KIAA1161 Polyclonal Antibody	A65678	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
KIAA1191 Polyclonal Antibody	A51141	Human, Mouse	ELISA, WB, IHC
KIAA1257 Polyclonal Antibody	A59550	Human	ELISA, IHC
KIAA1324L Polyclonal Antibody	A62822	Human	ELISA, IHC
KIAA1841 Polyclonal Antibody	A65510	Human	ELISA
KIF12 Polyclonal Antibody	A59554	Human	ELISA, IHC
KIF13B Polyclonal Antibody	A69231	Human	ELISA, IHC, IF
KIF15 Polyclonal Antibody	A69211	Human	ELISA, IF
Kif1a Polyclonal Antibody	A57085	Mouse	ELISA
KIF1BP Polyclonal Antibody	A70039	Human	ELISA, IHC, IF
Kif1c Polyclonal Antibody	A56663	Rat	ELISA
KIF20B Polyclonal Antibody	A69079	Human	ELISA, IHC, IF
KIF22 Polyclonal Antibody	A68859	Human	ELISA, IHC, IF
KIF24 Polyclonal Antibody	A65178	Human	ELISA, IHC
KIF26B Polyclonal Antibody	A64854	Human	ELISA, IHC
KIFAP3 Polyclonal Antibody	A68599	Human	ELISA, IF
KIFC1 Polyclonal Antibody	A70239	Human	ELISA, IF
KIR2DS1 Polyclonal Antibody	A57497	Human	ELISA
KIRREL2 Polyclonal Antibody	A65594	Human, Mouse	ELISA, WB
KLC4 Polyclonal Antibody	A62830	Human	ELISA
KLF1 Polyclonal Antibody	A64762	Human	ELISA
KLF13 Polyclonal Antibody	A70347	Human, Mouse, Rat	ELISA, WB, IHC
KLHDC8A Polyclonal Antibody	A59558	Human, Mouse	ELISA, WB, IHC
KLHL12 Polyclonal Antibody	A59562	Human	ELISA, IHC
KLHL2 Polyclonal Antibody	A62834	Human	ELISA
KLHL22 Polyclonal Antibody	A59566	Human, Mouse	ELISA, WB
KLHL23 Polyclonal Antibody	A66390	Human, Mouse	ELISA, WB
KLHL34 Polyclonal Antibody	A66942	Human	ELISA, IHC
KLHL38 Polyclonal Antibody	A64878	Human	ELISA, IHC
KLHL41 Polyclonal Antibody	A59570	Human	ELISA
KLHL42 Polyclonal Antibody	A59574	Human	ELISA, IHC, IF
KLK1 Polyclonal Antibody	A55243	Human	ELISA, IHC
KLK14 Polyclonal Antibody	A70231	Human	ELISA, IHC
Klk1b22 Polyclonal Antibody	A59578	Mouse	ELISA
KLK3 Polyclonal Antibody	A53035	Human	ELISA
KLK5 Polyclonal Antibody	A51429	Human	ELISA, WB, IHC
KLK6 Polyclonal Antibody	A66674	Human, Mouse	ELISA, WB, IHC, IF
Klk7 Polyclonal Antibody	A57121	Rat	ELISA
Klkb1 Polyclonal Antibody	A62838	Rat	ELISA
KLRC2 Polyclonal Antibody	A50326	Human	ELISA, IHC
KLRG1 Polyclonal Antibody	A55655	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
KMO Polyclonal Antibody	A55858	Human	ELISA, IHC, IF
KNG1 Polyclonal Antibody	A52162	Human	ELISA, WB, IHC, IF
Kng1 Polyclonal Antibody	A57109	Rat	ELISA
KP6 Killer Toxin Polyclonal Antibody	A64302	Ustilago maydis P6 virus	ELISA
KPNA1 Polyclonal Antibody	A54356	Human	ELISA, IHC
KRAS Polyclonal Antibody	A52273	Human, Mouse	ELISA, WB, IHC
KRBOX4 Polyclonal Antibody	A65158	Human, Mouse	ELISA, WB, IF
KRCC1 Polyclonal Antibody	A59582	Human, Mouse	ELISA, WB
KRT1 Polyclonal Antibody	A69567	Human	ELISA, IHC, IF
KRT10 Polyclonal Antibody	A54994	Human	ELISA, WB, IHC, IF
KRT14 Polyclonal Antibody	A56250	Human	ELISA, IHC, IF
KRT17 Polyclonal Antibody	A62842	Bovine	ELISA
KRT18 Polyclonal Antibody	A53182	Human, Zebrafish	ELISA, WB, IHC, IF
KRT19 Polyclonal Antibody	A52984	Human	ELISA
KRT19 Polyclonal Antibody	A60866	Human	ELISA, WB, IHC
KRT33B Polyclonal Antibody	A59586	Human	ELISA, WB, IHC
KRT35 Polyclonal Antibody	A67082	Human	ELISA, WB, IHC
KRT38 Polyclonal Antibody	A57141	Human	ELISA, WB
KRT4 Polyclonal Antibody	A55447	Human	ELISA, IHC, IF
KRT40 Polyclonal Antibody	A65190	Human	ELISA, IHC
Krt6a Polyclonal Antibody	A54535	Human	ELISA, IHC
KRT6A Polyclonal Antibody	A57041	Human	ELISA, WB, IHC, IF
KRT71 Polyclonal Antibody	A64926	Human	ELISA, IHC
KRT72 Polyclonal Antibody	A64774	Human	ELISA, WB
KRT73 Polyclonal Antibody	A65934	Human	ELISA, WB
KRT77 Polyclonal Antibody	A51009	Human	ELISA, IHC
KRT79 Polyclonal Antibody	A65214	Human, Mouse	ELISA, WB
KRT8 Polyclonal Antibody	A50676	Human, Zebrafish	ELISA, WB, IHC
KRT80 Polyclonal Antibody	A65642	Human, Mouse	ELISA, WB, IHC
KRT81 Polyclonal Antibody	A62846	Human, Mouse	ELISA, WB, IHC
KRTAP11-1 Polyclonal Antibody	A66058	Human	ELISA
KRTAP13-1 Polyclonal Antibody	A66134	Human	ELISA
KRTAP13-3 Polyclonal Antibody	A64954	Human	ELISA
KRTAP13-4 Polyclonal Antibody	A64958	Human	ELISA
KRTAP19-5 Polyclonal Antibody	A64910	Human	ELISA
KRTAP23-1 Polyclonal Antibody	A62850	Human	ELISA, IHC
KRTAP26-1 Polyclonal Antibody	A65698	Human	ELISA
KRTAP3-1 Polyclonal Antibody	A67234	Human	ELISA
KRTAP3-2 Polyclonal Antibody	A67458	Human	ELISA, IHC
KRTAP3-3 Polyclonal Antibody	A67550	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
KRTAP8-1 Polyclonal Antibody	A66114	Human	ELISA, IHC
KTN1 Polyclonal Antibody	A69971	Human, Mouse, Rat	ELISA, WB, IHC
KYAT1 Polyclonal Antibody	A56094	Human	ELISA, IHC
KYAT3 Polyclonal Antibody	A56086	Human	ELISA, IHC
L1 Polyclonal Antibody	A52150	Human papillomavirus type 18	ELISA, WB
L1TD1 Polyclonal Antibody	A65282	Human	ELISA, IHC
L2 Polyclonal Antibody	A56895	Human papillomavirus type 16	ELISA
L2 Polyclonal Antibody	A56907	Human papillomavirus type 18	ELISA
L2HGDH Polyclonal Antibody	A59590	Human	ELISA, WB, IHC
L2HGDH Polyclonal Antibody	A62854	Bovine	ELISA
L3HYPDH Polyclonal Antibody	A59594	Human, Mouse	ELISA, WB, IHC
L3MBTL4 Polyclonal Antibody	A66502	Human, Rat	ELISA, WB, IHC
LACE1 Polyclonal Antibody	A66258	Human, Mouse	ELISA, WB, IHC
LACTB2 Polyclonal Antibody	A59598	Human	ELISA, IHC, IF
Lactotransferrin Polyclonal Antibody	A57749	Human	ELISA
lacZ Polyclonal Antibody	A68539	Escherichia coli	ELISA
LAG3 Polyclonal Antibody	A56567	Human, Rat	ELISA, WB
LAIR1 Polyclonal Antibody	A62858	Human	ELISA, IHC, IF
LAIR2 Polyclonal Antibody	A62862	Human	ELISA, IHC
LAMA2 Polyclonal Antibody	A68603	Human	ELISA, IHC, IF
LAMA5 Polyclonal Antibody	A59602	Human	ELISA, IHC
LAMB2 Polyclonal Antibody	A52400	Human	ELISA
LAMC1 Polyclonal Antibody	A54628	Human	ELISA, IHC
LAMC2 Polyclonal Antibody	A68855	Human	ELISA, WB, IHC
LAMP2 Polyclonal Antibody	A50330	Human, Mouse	ELISA, WB, IHC, IF
LANCL2 Polyclonal Antibody	A70211	Human	ELISA, WB, IHC, IF
lapB Polyclonal Antibody	A53836	Escherichia coli	ELISA, WB
LAPTM4A Polyclonal Antibody	A69899	Human	ELISA, IHC, IF
LAPTM4B Polyclonal Antibody	A69987	Human	ELISA, WB, IHC, IF
LAPTM5 Polyclonal Antibody	A64814	Human	ELISA, IHC
LARP1B Polyclonal Antibody	A57133	Human	ELISA
LARP6 Polyclonal Antibody	A70319	Human	ELISA, IHC
LAS1L Polyclonal Antibody	A69351	Human	ELISA, WB, IF
lasA Polyclonal Antibody	A64266	Pseudomonas aeruginosa	ELISA
LCA5 Polyclonal Antibody	A65810	Human, Mouse	ELISA, WB
LCAT Polyclonal Antibody	A52082	Human	ELISA, WB
LCE2A Polyclonal Antibody	A65206	Human	ELISA, IHC
LCK Polyclonal Antibody	A56046	Human	ELISA, IHC, IF, IP

Product Name	Cat. #	Species Reactivity	Applications
LCN2 Polyclonal Antibody	A51525	Human	ELISA, WB
LCN2 Polyclonal Antibody	A51529	Human	ELISA, WB
LCN2 Polyclonal Antibody	A54029	Human	ELISA
Lcn3 Polyclonal Antibody	A65302	Mouse	ELISA
LCP1 Polyclonal Antibody	A50334	Human, Mouse	ELISA, WB, IHC, IF
LCP1 Polyclonal Antibody	A57157	Human	ELISA, WB
LCY1 Polyclonal Antibody	A57473	Arabidopsis thaliana	ELISA
LDAH Polyclonal Antibody	A59606	Human	ELISA, WB, IHC
IdcA Polyclonal Antibody	A53137	Escherichia coli	ELISA
IdcC Polyclonal Antibody	A52328	Escherichia coli	ELISA, WB
ldh1 Polyclonal Antibody	A54571	Lactobacillus plantarum	ELISA
LDHA Polyclonal Antibody	A52408	Human	ELISA, WB
LDHD Polyclonal Antibody	A65962	Human, Mouse	ELISA, WB, IHC, IF
LDLRAD1 Polyclonal Antibody	A65278	Human	ELISA, IF
lecA Polyclonal Antibody	A57189	Pseudomonas aeruginosa	ELISA
LECT1 Polyclonal Antibody	A59610	Human	ELISA
LEMD2 Polyclonal Antibody	A69143	Human	ELISA, WB, IHC
LEP Polyclonal Antibody	A53524	Human	ELISA, IHC, IF
lepA Polyclonal Antibody	A64418	Escherichia coli	ELISA
lepB Polyclonal Antibody	A64506	Escherichia coli	ELISA
lepB Polyclonal Antibody	A64602	Mycobacterium tuberculosis	ELISA
LGALS1 Polyclonal Antibody	A57097	Human	ELISA, WB, IHC, IF
LGALS12 Polyclonal Antibody	A59614	Human	ELISA, WB, IHC
LGALS14 Polyclonal Antibody	A66878	Human	ELISA
LGALS3 Polyclonal Antibody	A53624	Human	ELISA, WB, IHC
Lgals3bp Polyclonal Antibody	A54297	Rat	ELISA
LGALS7 Polyclonal Antibody	A52290	Human	ELISA
Lgals7 Polyclonal Antibody	A55786	Mouse	ELISA
LGALS9 Polyclonal Antibody	A55214	Human, Mouse	ELISA, WB, IHC
LGALS9C Polyclonal Antibody	A65266	Human, Mouse	ELISA, WB
LGI4 Polyclonal Antibody	A70023	Human	ELISA, IHC
LGMN Polyclonal Antibody	A69571	Human	ELISA, IHC
LHB Polyclonal Antibody	A54376	Human	ELISA, IHC
LHCGR Polyclonal Antibody	A50338	Human, Mouse	ELISA, WB, IHC, IF
Lhcgr Polyclonal Antibody	A50342	Rat	ELISA, WB
LHFP Polyclonal Antibody	A62866	Human	ELISA
LHFPL1 Polyclonal Antibody	A65922	Human	ELISA, IHC, IF
LHPP Polyclonal Antibody	A59618	Human	ELISA, IHC
LHX2 Polyclonal Antibody	A68607	Human, Rat	ELISA, WB, IF
LIFR Polyclonal Antibody	A69575	Human, Rat	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
LIG1 Polyclonal Antibody	A55399	Human	ELISA, IHC, IF
ligA Polyclonal Antibody	A53481	Escherichia coli	ELISA, WB
LILRA2 Polyclonal Antibody	A59622	Human	ELISA, WB, IHC
LILRA3 Polyclonal Antibody	A70059	Human, Mouse, Rat	ELISA, WB, IHC
LILRA4 Polyclonal Antibody	A59626	Human	ELISA, IHC, IF
LILRB1 Polyclonal Antibody	A70115	Human, Rat	ELISA, WB, IHC, IF
Lilrb3 Polyclonal Antibody	A55054	Mouse	ELISA
LIMS3 Polyclonal Antibody	A62870	Human	ELISA, IHC, IF
LIN37 Polyclonal Antibody	A59630	Human	ELISA, IHC
LIN7A Polyclonal Antibody	A62874	Human, Mouse, Zebrafish	ELISA, WB
LINC00846 Polyclonal Antibody	A59634	Human	ELISA
LINC01553 Polyclonal Antibody	A62162	Human, Mouse, Rat	ELISA, WB
LIPA Polyclonal Antibody	A53766	Human	ELISA, WB, IHC
LIPA Polyclonal Antibody	A54716	Human	ELISA, WB, IHC, IF
LIPH Polyclonal Antibody	A70127	Human	ELISA, IHC, IF
LIX1 Polyclonal Antibody	A59638	Human, Mouse	ELISA, WB, IHC
lktA Polyclonal Antibody	A57769	Mannheimia haemolytica	ELISA
L-Lactate Dehydrogenase Polyclonal Antibody	A55954	Plasmodium falciparum	ELISA
LLPH Polyclonal Antibody	A59642	Human	ELISA
LMAN1 Polyclonal Antibody	A68611	Human, Mouse	ELISA, WB, IHC, IF
LMAN1L Polyclonal Antibody	A56262	Human	ELISA, IHC
LMAN2L Polyclonal Antibody	A67246	Human	ELISA
LMBRD1 Polyclonal Antibody	A67862	Human, Mouse	ELISA, WB, IHC
LMCD1 Polyclonal Antibody	A70255	Human	ELISA, IHC, IF
LMF2 Polyclonal Antibody	A62878	Human, Mouse	ELISA, WB, IF
LMNTD1 Polyclonal Antibody	A59646	Human, Mouse	ELISA, WB
LMO1 Polyclonal Antibody	A53905	Human	ELISA, WB
LMO3 Polyclonal Antibody	A59650	Human	ELISA, WB, IHC
LNX2 Polyclonal Antibody	A67038	Human	ELISA, IHC
LOH12CR1 Polyclonal Antibody	A59654	Human	ELISA, IHC
lolA Polyclonal Antibody	A56899	Escherichia coli O9:H4	ELISA
LONRF1 Polyclonal Antibody	A64694	Human	ELISA, IHC
LOX Polyclonal Antibody	A52332	Human, Mouse	ELISA, IHC
Lox Polyclonal Antibody	A53578	Rat	ELISA, WB
Loxl1 Polyclonal Antibody	A52956	Human	ELISA, IHC
Loxl1 Polyclonal Antibody	A59658	Mouse	ELISA
Loxl2 Polyclonal Antibody	A56627	Mouse, Human	ELISA, WB
LOXL3 Polyclonal Antibody	A62882	Human	ELISA, IHC
LPA Polyclonal Antibody	A50833	Human	ELISA
LPAR4 Polyclonal Antibody	A62886	Human, Mouse	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
LPGAT1 Polyclonal Antibody	A66270	Human	ELISA, WB, IHC
lptA Polyclonal Antibody	A57237	Escherichia coli	ELISA
lpxC Polyclonal Antibody	A53321	Escherichia coli	ELISA
LPXN Polyclonal Antibody	A62890	Human	ELISA, WB, IHC, IF
LRBA Polyclonal Antibody	A52764	Human	ELISA, IHC, IF
LRG1 Polyclonal Antibody	A64136	Human, Mouse	ELISA, WB, IHC
LRP2 Polyclonal Antibody	A68615	Human	ELISA, IHC
LRPPRC Polyclonal Antibody	A69579	Human	ELISA, WB, IHC, IF
LRR15 Polyclonal Antibody	A66246	Human	ELISA, IHC
LRR18 Polyclonal Antibody	A66622	Human	ELISA, IHC
LRR19 Polyclonal Antibody	A67506	Human	ELISA, IHC
LRR2 Polyclonal Antibody	A59662	Human	ELISA, IHC
LRR20 Polyclonal Antibody	A66410	Human, Mouse	ELISA, WB, IHC
LRR29 Polyclonal Antibody	A66254	Human	ELISA
LRR31 Polyclonal Antibody	A65442	Human, Mouse	ELISA, WB, IHC
LRR32 Polyclonal Antibody	A64738	Human	ELISA, IHC
LRR39 Polyclonal Antibody	A59666	Human	ELISA, IHC
LRR3B Polyclonal Antibody	A67126	Human	ELISA, IHC, IF
LRR40 Polyclonal Antibody	A53788	Human	ELISA, WB
LRR42 Polyclonal Antibody	A68186	Human	ELISA, WB, IHC
LRR45 Polyclonal Antibody	A70131	Human	ELISA, IHC, IF
LRR4C Polyclonal Antibody	A67970	Human	ELISA, IHC, IF
LRR57 Polyclonal Antibody	A66762	Human, Mouse	ELISA, WB
LRR61 Polyclonal Antibody	A59670	Human	ELISA, IHC
LRR8E Polyclonal Antibody	A59674	Human	ELISA, IHC, IF
LRRFIP1 Polyclonal Antibody	A53344	Human	ELISA, IHC, IF
LRRN1 Polyclonal Antibody	A50997	Human	ELISA, IHC
LRRN2 Polyclonal Antibody	A62894	Human	ELISA, IHC, IF
LRRN3 Polyclonal Antibody	A67698	Human, Mouse	ELISA, WB, IHC, IF
LRRTM2 Polyclonal Antibody	A62898	Human	ELISA
LRTM1 Polyclonal Antibody	A67850	Human	ELISA, WB, IHC
LRTOMT Polyclonal Antibody	A52380	Human	ELISA, IF
LSAMP Polyclonal Antibody	A69583	Human, Mouse, Rat	ELISA, WB, IF
LSM1 Polyclonal Antibody	A62902	Human	ELISA, IHC
LSM10 Polyclonal Antibody	A55894	Human	ELISA, IHC
LSM3 Polyclonal Antibody	A59678	Human	ELISA
LSM4 Polyclonal Antibody	A50346	Human	ELISA, WB, IHC, IF
LSM6 Polyclonal Antibody	A59682	Human	ELISA
LSM7 Polyclonal Antibody	A68022	Human	ELISA, WB, IHC
LSMEM1 Polyclonal Antibody	A66202	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
LSMEM2 Polyclonal Antibody	A59686	Human	ELISA, IHC, IF
IspL Polyclonal Antibody	A64590	Rhizobium meliloti	ELISA
LSR Polyclonal Antibody	A69015	Human	ELISA, IF
LTB Polyclonal Antibody	A55543	Chimpanzee	ELISA
LTBR Polyclonal Antibody	A54077	Human	ELISA
LTF Polyclonal Antibody	A52456	Human	ELISA, WB
LTV1 Polyclonal Antibody	A67010	Human	ELISA, IHC
luxS Polyclonal Antibody	A54688	Escherichia coli	ELISA, WB
LUZP2 Polyclonal Antibody	A62906	Human	ELISA, WB, IHC, IF
LY6D Polyclonal Antibody	A64650	Human	ELISA
LY6E Polyclonal Antibody	A50917	Human	ELISA, IHC, IF
Ly6e Polyclonal Antibody	A59690	Mouse, Human	ELISA, WB
LY6G6C Polyclonal Antibody	A62910	Human	ELISA
LY6H Polyclonal Antibody	A62914	Human	ELISA, IHC, IF
Ly75 Polyclonal Antibody	A57437	Mouse	ELISA
LY96 Polyclonal Antibody	A59694	Human	ELISA, IHC
LYAR Polyclonal Antibody	A67750	Human	ELISA, WB, IHC
LYG1 Polyclonal Antibody	A59698	Human	ELISA, IHC
LYG2 Polyclonal Antibody	A65870	Human	ELISA, IHC
LYPD2 Polyclonal Antibody	A65654	Human	ELISA, IHC
LYPD3 Polyclonal Antibody	A62918	Human, Mouse	ELISA, WB
LYPD4 Polyclonal Antibody	A59702	Human	ELISA
Lypla1 Polyclonal Antibody	A53240	Mouse	ELISA
Lypla1 Polyclonal Antibody	A62922	Mouse, Human	ELISA, IF
LYPLA2 Polyclonal Antibody	A53989	Human	ELISA, IF
LYPLAL1 Polyclonal Antibody	A59706	Human	ELISA, IHC, IF
LYRM1 Polyclonal Antibody	A59710	Human	ELISA, IHC
LYRM2 Polyclonal Antibody	A67858	Human	ELISA, IHC
LYRM5 Polyclonal Antibody	A65558	Human	ELISA, IHC
LYSMD1 Polyclonal Antibody	A66474	Human	ELISA, IHC
LYSMD3 Polyclonal Antibody	A65786	Human, Mouse	ELISA, WB
LYZ Polyclonal Antibody	A50616	Human	ELISA, IHC
LYZL1 Polyclonal Antibody	A50993	Human	ELISA, IHC
LYZL2 Polyclonal Antibody	A65894	Human	ELISA, IHC
LZTS1 Polyclonal Antibody	A68783	Human	ELISA, IHC, IF
M6PR Polyclonal Antibody	A53316	Human	ELISA, IHC, IF
MAB21L1 Polyclonal Antibody	A64838	Human, Mouse	ELISA, WB, IHC
MACROD1 Polyclonal Antibody	A59714	Human	ELISA, IHC
MAD2L1 Polyclonal Antibody	A50921	Human	ELISA, IHC
MAD2L2 Polyclonal Antibody	A59718	Human	ELISA



Product Name	Cat. #	Species Reactivity	Applications
MADCAM1 Polyclonal Antibody	A64726	Human	ELISA, WB, IF
MAEL Polyclonal Antibody	A68266	Human	ELISA, IHC
MAF1 Polyclonal Antibody	A56074	Human, Mouse	ELISA, WB
MAG Polyclonal Antibody	A54412	Human	ELISA
MAGEA10 Polyclonal Antibody	A50350	Human	ELISA, WB, IHC
MAGEA4 Polyclonal Antibody	A59726	Human	ELISA, IHC
MAGEA5 Polyclonal Antibody	A62926	Human	ELISA, IHC
MAGEA8 Polyclonal Antibody	A50354	Human	ELISA, IHC, IF
MAGEA9 Polyclonal Antibody	A59730	Human	ELISA, IHC
MAGEB10 Polyclonal Antibody	A67026	Human, Mouse	ELISA, WB
MAGEB18 Polyclonal Antibody	A59734	Human, Mouse	ELISA, WB
MAGEC3 Polyclonal Antibody	A59738	Human, Rat	ELISA, WB
MAGED2 Polyclonal Antibody	A68090	Human	ELISA
MAGOHB Polyclonal Antibody	A59742	Human	ELISA, IHC
Major Surface Glycoprotein G Polyclonal Antibody	A50853	Human respiratory syncytial virus A	ELISA
MAMDC2 Polyclonal Antibody	A59746	Human	ELISA
manA Polyclonal Antibody	A54998	Escherichia coli	ELISA, WB
MANBAL Polyclonal Antibody	A59750	Human	ELISA, IHC, IF
MAOA Polyclonal Antibody	A59754	Human, Mouse	ELISA, WB, IHC
MAOB Polyclonal Antibody	A53194	Human, Mouse	ELISA, WB, IHC
MAP10 Polyclonal Antibody	A67646	Human, Mouse	ELISA, WB, IHC, IF
MAP1A Polyclonal Antibody	A68619	Human	ELISA, IHC, IF
MAP1LC3A Polyclonal Antibody	A51349	Human	ELISA, IHC
MAP1S Polyclonal Antibody	A52253	Human	ELISA, IHC
MAP2K3 Polyclonal Antibody	A54584	Human, Mouse	ELISA, WB, IHC, IF
Map2k4 Polyclonal Antibody	A56503	Mouse, Human	ELISA, WB
MAP2K6 Polyclonal Antibody	A50358	Human	ELISA, WB
MAP3K1 Polyclonal Antibody	A52118	Mouse	ELISA
MAP3K12 Polyclonal Antibody	A62930	Human	ELISA
MAP3K19 Polyclonal Antibody	A65118	Human	ELISA
MAP3K7 Polyclonal Antibody	A62934	Human	ELISA, IHC
MAP3K7CL Polyclonal Antibody	A50809	Human	ELISA, IHC
MAP3K8 Polyclonal Antibody	A62938	Human	ELISA, IHC, IF
MAPK13 Polyclonal Antibody	A59758	Human	ELISA, WB, IHC, IF
MAPK3 Polyclonal Antibody	A54955	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC, IF
MAPK9 Polyclonal Antibody	A50362	Human, Mouse	ELISA, WB, IP
MAPKAPK5 Polyclonal Antibody	A66158	Human	ELISA
MAPRE1 Polyclonal Antibody	A64782	Human	ELISA, IHC
Mapt Polyclonal Antibody	A54812	Rat	ELISA

Product Name	Cat. #	Species Reactivity	Applications
MAS1 Polyclonal Antibody	A50366	Human, Mouse	ELISA, WB, IHC, IF
MAS1L Polyclonal Antibody	A59762	Human	ELISA, WB, IHC, IF
MASP1 Polyclonal Antibody	A59766	Human	ELISA, WB
MAT2A Polyclonal Antibody	A62942	Human	ELISA, IHC, IF
Matrix protein Polyclonal Antibody	A54616	Rabies virus	ELISA, WB
MB Polyclonal Antibody	A54065	Human, Mouse, Rat	ELISA, WB, IHC
MB Polyclonal Antibody	A54256	Human, Mouse, Rat	ELISA, WB, IHC
MBD5 Polyclonal Antibody	A67422	Human	ELISA
MBIP Polyclonal Antibody	A70323	Human	ELISA, IHC
MBNL1 Polyclonal Antibody	A55535	Human	ELISA, IHC
MC3R Polyclonal Antibody	A69587	Human	ELISA, IHC
MCAM Polyclonal Antibody	A53105	Human	ELISA, IHC
Mcl1 Polyclonal Antibody	A55094	Mouse	ELISA
MCM2 Polyclonal Antibody	A54744	Human	ELISA, WB
MCM9 Polyclonal Antibody	A59770	Human, Mouse	ELISA, WB, IHC
Mcpt4 Polyclonal Antibody	A54890	Mouse	ELISA
MCTS1 Polyclonal Antibody	A51401	Human	ELISA, IHC
MDK Polyclonal Antibody	A52856	Human, Mouse	ELISA, IHC, IF
MDR1 Polyclonal Antibody	A50740	Plasmodium falciparum	ELISA
MDR1 Polyclonal Antibody	A62034	Entamoeba histolytica	ELISA
MECR Polyclonal Antibody	A51369	Human, Rat	ELISA, WB, IHC
MED15 Polyclonal Antibody	A69091	Human	ELISA, WB, IF
MED18 Polyclonal Antibody	A59774	Human	ELISA, WB, IHC
MED22 Polyclonal Antibody	A64778	Human	ELISA, IHC, IF
MED24 Polyclonal Antibody	A62946	Human	ELISA, IHC
MED27 Polyclonal Antibody	A65690	Human	ELISA, WB
MED28 Polyclonal Antibody	A59778	Human	ELISA, IHC
MED30 Polyclonal Antibody	A59782	Human	ELISA, IHC, IF
MED7 Polyclonal Antibody	A59786	Human, Mouse	ELISA, WB
MEGF9 Polyclonal Antibody	A69223	Human	ELISA, WB, IHC, IF
MEI1 Polyclonal Antibody	A59790	Human	ELISA, IHC
MEIS3 Polyclonal Antibody	A68623	Human	ELISA, IHC
meIA Polyclonal Antibody	A54712	Bacillus subtilis	ELISA
MELTF Polyclonal Antibody	A62954	Human	ELISA, WB, IHC, IF
METAP1 Polyclonal Antibody	A68627	Human, Rat	ELISA, WB, IHC, IF
metC Polyclonal Antibody	A53405	Escherichia coli	ELISA
METRNL Polyclonal Antibody	A56114	Human, Mouse, Rat	ELISA, WB, IHC
METTL25 Polyclonal Antibody	A66630	Human	ELISA, IHC
METTL2A Polyclonal Antibody	A67102	Human, Mouse	ELISA, WB
METTL7A Polyclonal Antibody	A67198	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
MFAP3 Polyclonal Antibody	A62950	Human	ELISA
MFAP3L Polyclonal Antibody	A59794	Human	ELISA, IHC, IF
MFAP4 Polyclonal Antibody	A69591	Human	ELISA, IHC
MFAP5 Polyclonal Antibody	A56134	Human	ELISA, IHC
MFGE8 Polyclonal Antibody	A53057	Human, Mouse	ELISA, WB, IHC
MFSD1 Polyclonal Antibody	A67262	Human	ELISA, IHC, IF
MFSD11 Polyclonal Antibody	A62958	Human	ELISA, IHC
MFSD14B Polyclonal Antibody	A59370	Human	ELISA, IHC, IF
MFSD5 Polyclonal Antibody	A59798	Human	ELISA, IHC, IF
MFSD8 Polyclonal Antibody	A66782	Human	ELISA, IHC, IF
MGAT3 Polyclonal Antibody	A69799	Human, Mouse	ELISA, WB, IHC
MGME1 Polyclonal Antibody	A59802	Human	ELISA, WB, IHC
MGMT Polyclonal Antibody	A52416	Human	ELISA, IHC
MGP Polyclonal Antibody	A57417	Human	ELISA
mgsA Polyclonal Antibody	A56238	Bacillus cereus	ELISA
MHT1 Polyclonal Antibody	A52209	Saccharomyces cerevisiae	ELISA, WB
MIA3 Polyclonal Antibody	A65150	Human	ELISA
MIB1 Polyclonal Antibody	A54177	Human	ELISA, WB, IHC, IF
MICAL2 Polyclonal Antibody	A59806	Human	ELISA, IHC, IF
MICU1 Polyclonal Antibody	A70179	Human	ELISA, WB, IHC
MIEF1 Polyclonal Antibody	A59810	Human, Mouse	ELISA, WB
MIER2 Polyclonal Antibody	A59814	Human	ELISA, WB, IHC
MIF Polyclonal Antibody	A51917	Human	ELISA, WB
MIF4GD Polyclonal Antibody	A59818	Human	ELISA
Milk Beta-Lactoglobulin Polyclonal Antibody	A57781	Bovine	ELISA, WB
MINA Polyclonal Antibody	A66062	Human	ELISA, IHC, IF
Minor Capsid Protein VP2 Polyclonal Antibody	A55206	JC polyomavirus	ELISA
Minpp1 Polyclonal Antibody	A50370	Rat	ELISA
Minpp1 Polyclonal Antibody	A50374	Rat	ELISA
MINPP1 Polyclonal Antibody	A56126	Human	ELISA, WB, IHC
MIOX Polyclonal Antibody	A67338	Human, Mouse	ELISA, WB
mip Polyclonal Antibody	A64562	Legionella pneumophila	ELISA
MIP Polyclonal Antibody	A69595	Human	ELISA, IHC, IF
MIS12 Polyclonal Antibody	A51329	Human	ELISA, WB, IHC
MKI67 Polyclonal Antibody	A53501	Human	ELISA, IHC
MKX Polyclonal Antibody	A70019	Human	ELISA, WB, IHC, IF
MLANA Polyclonal Antibody	A50929	Human	ELISA, WB, IHC
MLKL Polyclonal Antibody	A51161	Human	ELISA, IHC
MLLT10 Polyclonal Antibody	A62962	Human	ELISA, IHC
MLLT11 Polyclonal Antibody	A59822	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
MLX Polyclonal Antibody	A59826	Human	ELISA, IHC, IF
MLXIP Polyclonal Antibody	A51237	Human	ELISA, IHC
MMADHC Polyclonal Antibody	A70267	Human	ELISA, IHC, IF
MMD2 Polyclonal Antibody	A66126	Human, Mouse	ELISA, WB, IHC, IF
MMGT1 Polyclonal Antibody	A50378	Human	ELISA, WB
MMP1 Polyclonal Antibody	A54041	Human	ELISA, WB, IHC
MMP10 Polyclonal Antibody	A53300	Human, Mouse	ELISA, WB, IHC, IF
MMP12 Polyclonal Antibody	A54464	Human	ELISA, WB, IHC
Mmp13 Polyclonal Antibody	A52508	Mouse, Human	ELISA, IHC
MMP13 Polyclonal Antibody	A55291	Human, Mouse	ELISA, WB, IHC, IF
MMP14 Polyclonal Antibody	A55030	Human	ELISA
Mmp2 Polyclonal Antibody	A57113	Rat	ELISA
MMP2 Polyclonal Antibody	A64096	Human	ELISA, IHC, IF
Mmp3 Polyclonal Antibody	A52512	Mouse, Human	ELISA, WB, IHC
MMP3 Polyclonal Antibody	A55223	Human, Mouse	ELISA, WB, IHC
MMP7 Polyclonal Antibody	A51829	Human	ELISA, IHC
Mmp7 Polyclonal Antibody	A53820	Rat, Mouse, Human	ELISA, WB
MND1 Polyclonal Antibody	A67662	Human	ELISA, WB, IHC
MOB1A Polyclonal Antibody	A57065	Human, Mouse	ELISA, WB, IHC
MOB1B Polyclonal Antibody	A56110	Human	ELISA, WB
MOBP Polyclonal Antibody	A69599	Human	ELISA, WB, IF
MOCS2 Polyclonal Antibody	A62966	Human	ELISA, WB
MOGAT2 Polyclonal Antibody	A64946	Human	ELISA, IHC
MON1A Polyclonal Antibody	A56034	Human	ELISA, WB, IHC
MORC2 Polyclonal Antibody	A59830	Human	ELISA, WB, IHC
MORN3 Polyclonal Antibody	A65482	Human, Mouse	ELISA, WB
MORN5 Polyclonal Antibody	A65358	Human, Mouse	ELISA, WB, IHC
Mouse Hemoglobin Polyclonal Antibody	A57573	Mouse	ELISA, WB
Mouse Serum Albumin Polyclonal Antibody	A57545	Mouse	ELISA
MOV10 Polyclonal Antibody	A68294	Human	ELISA, WB, IHC
MP68 Polyclonal Antibody	A51636	Human	ELISA, WB, IHC, IF
MPDU1 Polyclonal Antibody	A62970	Human	ELISA, IHC, IF
MPO Polyclonal Antibody	A54284	Human	ELISA, IF
MPPED1 Polyclonal Antibody	A59834	Human	ELISA, IHC, IF
MPPED2 Polyclonal Antibody	A59838	Human, Mouse	ELISA, WB
mpt63 Polyclonal Antibody	A55363	Mycobacterium tuberculosis	ELISA, WB
mpt63 Polyclonal Antibody	A57627	Mycobacterium tuberculosis	ELISA
mpt64 Polyclonal Antibody	A55719	Mycobacterium tuberculosis	ELISA
mpt64 Polyclonal Antibody	A57629	Mycobacterium tuberculosis	ELISA, WB
MPV17 Polyclonal Antibody	A57269	Human, Mouse	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
MPZ Polyclonal Antibody	A52676	Human	ELISA, WB
MPZ Polyclonal Antibody	A54061	Human	ELISA
MPZL2 Polyclonal Antibody	A62974	Human	ELISA, IHC
MRAP2 Polyclonal Antibody	A67006	Human, Mouse	ELISA, WB, IHC, IF
MRC1 Polyclonal Antibody	A59842	Human	ELISA
MREG Polyclonal Antibody	A62978	Human, Mouse	ELISA, WB
MRFAP1 Polyclonal Antibody	A68198	Human	ELISA
MRFAP1L1 Polyclonal Antibody	A53332	Human	ELISA, WB
MRGPRX2 Polyclonal Antibody	A66606	Human, Mouse	ELISA, WB, IF
MRGPRX3 Polyclonal Antibody	A66926	Human	ELISA, IF
MRJP1 Polyclonal Antibody	A53730	Apis mellifera	ELISA
MRM2 Polyclonal Antibody	A67782	Human	ELISA, IHC
MRM3 Polyclonal Antibody	A60662	Human	ELISA, IHC
MRPL1 Polyclonal Antibody	A56871	Human	ELISA, WB, IHC
MRPL11 Polyclonal Antibody	A68170	Human	ELISA, WB, IF
MRPL14 Polyclonal Antibody	A62982	Human, Mouse	ELISA, WB, IHC
MRPL16 Polyclonal Antibody	A59846	Human	ELISA, IHC
MRPL18 Polyclonal Antibody	A59850	Human	ELISA, WB, IHC
MRPL19 Polyclonal Antibody	A53781	Human, Mouse	ELISA, WB, IHC, IF
MRPL2 Polyclonal Antibody	A59854	Human	ELISA, IHC
MRPL20 Polyclonal Antibody	A55878	Human	ELISA, IHC
MRPL21 Polyclonal Antibody	A65890	Human	ELISA, IHC
MRPL22 Polyclonal Antibody	A59858	Human	ELISA, WB, IHC
MRPL27 Polyclonal Antibody	A67418	Human	ELISA, IHC
MRPL3 Polyclonal Antibody	A59862	Human	ELISA, WB, IHC
MRPL30 Polyclonal Antibody	A66238	Human	ELISA
MRPL33 Polyclonal Antibody	A59866	Human	ELISA, IHC
MRPL36 Polyclonal Antibody	A67874	Human	ELISA
MRPL38 Polyclonal Antibody	A59870	Human	ELISA, IHC
MRPL41 Polyclonal Antibody	A66086	Human	ELISA, IHC
MRPL42 Polyclonal Antibody	A55886	Human	ELISA
MRPL44 Polyclonal Antibody	A59874	Human	ELISA, WB, IHC
MRPL46 Polyclonal Antibody	A67694	Human	ELISA
MRPL48 Polyclonal Antibody	A59878	Human, Mouse	ELISA, WB, IHC
MRPL49 Polyclonal Antibody	A64638	Human, Mouse	ELISA, WB, IHC
MRPL50 Polyclonal Antibody	A59882	Human	ELISA, IHC
MRPL51 Polyclonal Antibody	A65066	Human	ELISA, WB, IHC
MRPL53 Polyclonal Antibody	A59886	Human	ELISA, IHC
MRPL54 Polyclonal Antibody	A54692	Human	ELISA, IHC
MRPL55 Polyclonal Antibody	A53785	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
MRPL9 Polyclonal Antibody	A51837	Human	ELISA, WB
MRPS10 Polyclonal Antibody	A59890	Human	ELISA, IHC
MRPS14 Polyclonal Antibody	A59894	Human	ELISA
MRPS18A Polyclonal Antibody	A67742	Human	ELISA, WB
MRPS18C Polyclonal Antibody	A59898	Human	ELISA
MRPS2 Polyclonal Antibody	A59902	Human	ELISA, WB
MRPS28 Polyclonal Antibody	A59906	Human	ELISA, IHC
MRPS31 Polyclonal Antibody	A59910	Human	ELISA, WB, IHC
MRPS36 Polyclonal Antibody	A59914	Human	ELISA, IHC
MRPS5 Polyclonal Antibody	A59918	Human	ELISA, WB
MRRF Polyclonal Antibody	A59922	Human	ELISA, IHC
Ms4a1 Polyclonal Antibody	A54836	Mouse, Human	ELISA, WB
MS4A3 Polyclonal Antibody	A59926	Human	ELISA, IF
MS4A4A Polyclonal Antibody	A66310	Human	ELISA, IHC, IF
MS4A6E Polyclonal Antibody	A67090	Human	ELISA, IHC
MS4A7 Polyclonal Antibody	A59930	Human	ELISA, IHC, IF
MS4A8 Polyclonal Antibody	A67798	Human	ELISA, IHC, IF
MSA2 PFB0300c Polyclonal Antibody	A57709	Plasmodium falciparum	ELISA
MSH4 Polyclonal Antibody	A62986	Human, Mouse	ELISA, WB, IHC, IF
MSH6 Polyclonal Antibody	A50386	Human	ELISA, IHC, IF
Msln Polyclonal Antibody	A55555	Mouse	ELISA
MSLN Polyclonal Antibody	A55563	Human	ELISA, IHC
MSTN Polyclonal Antibody	A54157	Human, Mouse	ELISA, WB, IHC
MSTO1 Polyclonal Antibody	A59934	Human	ELISA, IHC
MT1A Polyclonal Antibody	A57677	Human	ELISA, IHC
MT1B Polyclonal Antibody	A64188	Human	ELISA
MT1E Polyclonal Antibody	A57577	Human	ELISA, IHC, IF
MT1F Polyclonal Antibody	A64184	Human	ELISA
MT1G Polyclonal Antibody	A57673	Human	ELISA, IHC
MT1H Polyclonal Antibody	A59938	Human	ELISA
MT1M Polyclonal Antibody	A59942	Human	ELISA
MT1X Polyclonal Antibody	A57661	Human	ELISA, IHC
MT2731 Polyclonal Antibody	A64200	Mycobacterium tuberculosis	ELISA
MT2A Polyclonal Antibody	A57676	Human	ELISA, IHC
MT3 Polyclonal Antibody	A68755	Human	ELISA, IF
MT4 Polyclonal Antibody	A64192	Human	ELISA
MTBP Polyclonal Antibody	A70171	Human	ELISA, IHC
MTHFD2L Polyclonal Antibody	A59946	Human, Mouse	ELISA, WB
MTHFR Polyclonal Antibody	A50390	Human	ELISA, WB, IHC, IF
MTHFR Polyclonal Antibody	A59950	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
MTHFS Polyclonal Antibody	A50394	Human	ELISA, WB, IHC, IF
MTMR11 Polyclonal Antibody	A62990	Human	ELISA
MTMR6 Polyclonal Antibody	A59954	Human	ELISA
MTMR7 Polyclonal Antibody	A56531	Human, Mouse	ELISA, WB
MTR Polyclonal Antibody	A59958	Human	ELISA, IHC
MTRR Polyclonal Antibody	A51385	Human	ELISA, WB, IHC
MUC1 Polyclonal Antibody	A52829	Human	ELISA, IHC, IF
MUC16 Polyclonal Antibody	A54125	Human	ELISA, IHC
MUC2 Polyclonal Antibody	A62994	Human	ELISA
MUC20 Polyclonal Antibody	A69063	Human	ELISA, WB, IF
MUCL1 Polyclonal Antibody	A67814	Human	ELISA, IHC
Mug1 Polyclonal Antibody	A57389	Mouse	ELISA
Mup11 Polyclonal Antibody	A54563	Mouse	ELISA, WB
Mup11 Polyclonal Antibody	A54804	Mouse	ELISA
Mup2 Polyclonal Antibody	A54800	Mouse	ELISA
Mup2 Polyclonal Antibody	A54820	Mouse	ELISA
MUS81 Polyclonal Antibody	A70047	Human	ELISA, IHC, IF
mutT Polyclonal Antibody	A64466	Escherichia coli	ELISA
MVD Polyclonal Antibody	A69603	Human, Rat	ELISA, WB, IHC, IF
MVP Polyclonal Antibody	A56483	Human, Rat	ELISA, WB
MX2 Polyclonal Antibody	A69607	Human	ELISA, IHC, IF
MXD3 Polyclonal Antibody	A59962	Human	ELISA, IHC
MYADM Polyclonal Antibody	A59966	Human	ELISA, IHC, IF
Mybpc3 Polyclonal Antibody	A54926	Rat	ELISA
Myc Polyclonal Antibody	A53590	Mouse	ELISA
MYC Polyclonal Antibody	A63788	Human, Mouse	ELISA, WB, IHC
MYCBP Polyclonal Antibody	A51221	Human	ELISA, IHC
MYEF2 Polyclonal Antibody	A55938	Human	ELISA, IHC, IF
MYH11 Polyclonal Antibody	A69611	Human	ELISA, IHC
MYH4 Polyclonal Antibody	A68791	Human	ELISA, IHC, IF
MYH7B Polyclonal Antibody	A59970	Human	ELISA, IHC
MYH9 Polyclonal Antibody	A53889	Human	ELISA, IHC, IF
MYL2 Polyclonal Antibody	A64120	Human	ELISA, IHC, IF
MYL6 Polyclonal Antibody	A53150	Human	ELISA, IHC, IF
MYLPF Polyclonal Antibody	A56827	Human, Mouse	ELISA, WB
MYNN Polyclonal Antibody	A59974	Human	ELISA, IHC
MYO19 Polyclonal Antibody	A51081	Human	ELISA, WB
N4BP2L2 Polyclonal Antibody	A66570	Human	ELISA, IHC
N6AMT2 Polyclonal Antibody	A51061	Human	ELISA, WB
NAA38 Polyclonal Antibody	A67434	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
NAA60 Polyclonal Antibody	A70271	Human	ELISA, IHC
NAALADL2 Polyclonal Antibody	A56611	Human, Mouse	ELISA, WB, IHC, IF
NABP1 Polyclonal Antibody	A55838	Human	ELISA, WB
NACA2 Polyclonal Antibody	A67934	Human	ELISA, WB
NADK Polyclonal Antibody	A63002	Human	ELISA, WB
NADK2 Polyclonal Antibody	A69919	Human	ELISA, WB, IHC
NAGK Polyclonal Antibody	A56735	Human	ELISA, WB
NAIF1 Polyclonal Antibody	A59978	Human, Mouse	ELISA, WB, IHC
NANOG Polyclonal Antibody	A59982	Human	ELISA, WB, IHC
NANOGP8 Polyclonal Antibody	A65570	Human, Mouse, Rat	ELISA, WB
NANOS2 Polyclonal Antibody	A63006	Human	ELISA, IHC, IF
NANP Polyclonal Antibody	A67054	Human	ELISA, IHC
NANS Polyclonal Antibody	A69287	Human	ELISA, WB, IF
NAPA Polyclonal Antibody	A59986	Human	ELISA, WB, IHC, IF
NAPEPLD Polyclonal Antibody	A65674	Human, Mouse	ELISA, WB
NAPRT Polyclonal Antibody	A53726	Human, Mouse	ELISA, WB
NAPSA Polyclonal Antibody	A54491	Human, Mouse	ELISA, WB, IHC
NARFL Polyclonal Antibody	A67278	Human	ELISA, WB, IHC, IF
NARS Polyclonal Antibody	A63010	Human	ELISA, IHC, IF
NARS2 Polyclonal Antibody	A59990	Human	ELISA, IHC
NAT2 Polyclonal Antibody	A59994	Human, Mouse	ELISA, WB, IHC, IF
NAT8 Polyclonal Antibody	A59998	Human	ELISA, IHC, IF
NBPF3 Polyclonal Antibody	A67478	Human	ELISA, WB, IHC
NCDN Polyclonal Antibody	A69615	Human, Rat	ELISA, WB, IHC, IF
NCK2 Polyclonal Antibody	A56058	Human, Mouse	ELISA, WB, IHC, IF
NCMAP Polyclonal Antibody	A65346	Human	ELISA, IHC, IF
NCOR1 Polyclonal Antibody	A69619	Human	ELISA, IF
ND-75 Polyclonal Antibody	A66278	Drosophila melanogaster	ELISA
NDFIP1 Polyclonal Antibody	A60002	Human	ELISA, WB, IHC, IF
NDN Polyclonal Antibody	A70175	Human	ELISA, IHC, IF
NDST1 Polyclonal Antibody	A55978	Human	ELISA, WB, IHC
NDUFA1 Polyclonal Antibody	A51437	Human	ELISA, IHC, IF
NDUFA2 Polyclonal Antibody	A53089	Human	ELISA, IHC
NDUFA3 Polyclonal Antibody	A54208	Human	ELISA, IHC, IF
NDUFA6 Polyclonal Antibody	A54232	Human	ELISA, IHC
NDUFAB1 Polyclonal Antibody	A68210	Human	ELISA, IHC
NDUFAB2 Polyclonal Antibody	A69071	Human	ELISA, WB, IHC
NDUFB11 Polyclonal Antibody	A60006	Human	ELISA, IHC, IF
NDUFB5 Polyclonal Antibody	A50398	Human, Mouse	ELISA, WB
NDUFB5 Polyclonal Antibody	A60242	Human	ELISA



Product Name	Cat. #	Species Reactivity	Applications
NDUFB6 Polyclonal Antibody	A63014	Human, Mouse	ELISA, WB, IHC, IP
NDUFB7 Polyclonal Antibody	A52022	Human	ELISA, IHC
NDUFB8 Polyclonal Antibody	A69623	Human, Mouse, Rat	ELISA, WB, IHC
NDUFS5 Polyclonal Antibody	A51632	Human	ELISA, WB
NDUFV2 Polyclonal Antibody	A51661	Human	ELISA, WB
NDUFV3 Polyclonal Antibody	A63018	Human	ELISA
NEB Polyclonal Antibody	A69819	Human	ELISA, IHC, IF
NEGR1 Polyclonal Antibody	A60010	Human	ELISA, WB, IHC, IF
NEIL1 Polyclonal Antibody	A55739	Human	ELISA, IHC
NEK1 Polyclonal Antibody	A70051	Human	ELISA, IHC, IF
NEK11 Polyclonal Antibody	A60014	Human	ELISA, WB, IHC
NEK4 Polyclonal Antibody	A68631	Human	ELISA, IF
NEK5 Polyclonal Antibody	A69927	Human	ELISA, IHC
NELFB Polyclonal Antibody	A68308	Human	ELISA, WB, IHC, ChIP
NET1 Polyclonal Antibody	A55443	Human	ELISA
NETO1 Polyclonal Antibody	A70083	Human	ELISA, WB, IF
NETO2 Polyclonal Antibody	A66214	Human	ELISA, WB, IHC
NEU1 Polyclonal Antibody	A64180	Human	ELISA, IHC, IF
NEU4 Polyclonal Antibody	A70123	Human	ELISA, IHC, IF
NFAM1 Polyclonal Antibody	A66974	Human	ELISA, IHC
NFE2L1 Polyclonal Antibody	A63022	Human	ELISA, IHC, IF
NFIC Polyclonal Antibody	A63026	Human	ELISA, IHC, IF
NFIX Polyclonal Antibody	A64710	Human	ELISA, IHC
NFKB1 Polyclonal Antibody	A68309	Human	ELISA, WB, IHC, ChIP
NFKBID Polyclonal Antibody	A66786	Human	ELISA, WB
NGB Polyclonal Antibody	A55455	Human	ELISA
Ngf Polyclonal Antibody	A52042	Rat	ELISA
NGF Polyclonal Antibody	A64862	Pig	ELISA
NHLRC2 Polyclonal Antibody	A60018	Human	ELISA, WB, IHC
NHP2 Polyclonal Antibody	A67522	Human	ELISA, IHC, IF
NICN1 Polyclonal Antibody	A67342	Human	ELISA, WB, IHC
Nid1 Polyclonal Antibody	A54531	Mouse, Rat	ELISA, WB
Nid1 Polyclonal Antibody	A55070	Mouse	ELISA
NID1 Polyclonal Antibody	A68635	Human	ELISA, IHC
NIF3L1 Polyclonal Antibody	A60022	Human	ELISA, WB, IHC
nikR Polyclonal Antibody	A57001	Escherichia coli	ELISA
NIM1K Polyclonal Antibody	A54360	Human	ELISA, WB, IHC
NINJ1 Polyclonal Antibody	A67086	Human	ELISA, IHC, IF
NINJ2 Polyclonal Antibody	A67870	Human	ELISA, IHC, IF
NIP7 Polyclonal Antibody	A51425	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
NIPSNAP3B Polyclonal Antibody	A67658	Human	ELISA, IHC
NIT1 Polyclonal Antibody	A65766	Human	ELISA, IHC
NIT2 Polyclonal Antibody	A67726	Human, Mouse	ELISA, WB
NKAIN4 Polyclonal Antibody	A66154	Human	ELISA, IHC, IF
NKD1 Polyclonal Antibody	A70151	Human, Rat	ELISA, WB, IF
NKG7 Polyclonal Antibody	A60026	Human	ELISA, IF
NKIRAS1 Polyclonal Antibody	A60030	Human	ELISA, WB, IHC
NKIRAS2 Polyclonal Antibody	A60034	Human	ELISA, IHC
NKTR Polyclonal Antibody	A54320	Human	ELISA
Nkx2-2 Polyclonal Antibody	A60038	Mouse	ELISA
NKX2-8 Polyclonal Antibody	A69627	Human	ELISA, IHC
NLE1 Polyclonal Antibody	A51381	Human	ELISA
NLGN4X Polyclonal Antibody	A69099	Human	ELISA, IF
NLRC5 Polyclonal Antibody	A69967	Human	ELISA, IHC, IF
NLRP1 Polyclonal Antibody	A67466	Human	ELISA
NLRP10 Polyclonal Antibody	A68995	Human, Rat	ELISA, WB, IHC, IF
NLRP7 Polyclonal Antibody	A69087	Human	ELISA, IHC
NME1 Polyclonal Antibody	A53929	Human	ELISA, WB, IHC
NME5 Polyclonal Antibody	A63030	Human	ELISA, IF
NME6 Polyclonal Antibody	A50402	Human	ELISA, WB, IHC
NME7 Polyclonal Antibody	A60042	Human, Mouse	ELISA, WB, IHC
NMNAT3 Polyclonal Antibody	A57005	Human, Mouse	ELISA, WB, IHC
NMRAL1 Polyclonal Antibody	A60046	Human	ELISA, WB
NMUR1 Polyclonal Antibody	A67838	Human, Mouse	ELISA, WB, IHC, IF
NOL10 Polyclonal Antibody	A60050	Human	ELISA, WB, IHC
NOL3 Polyclonal Antibody	A56839	Human, Mouse	ELISA, WB
NOL4L Polyclonal Antibody	A67118	Human	ELISA, WB, IHC, IF
NOL6 Polyclonal Antibody	A60054	Human	ELISA, IHC
NOMO1 Polyclonal Antibody	A69631	Human	ELISA, WB, IHC, IF
NOP10 Polyclonal Antibody	A67518	Human	ELISA, WB, IHC
NOP58 Polyclonal Antibody	A54768	Human, Mouse	ELISA, WB, IHC, IF
Nos3 Polyclonal Antibody	A52138	Rat	ELISA
NOSTRIN Polyclonal Antibody	A60058	Human, Mouse	ELISA, WB, IHC
Nox4 Polyclonal Antibody	A54539	Rat	ELISA
NOX4 Polyclonal Antibody	A63034	Human	ELISA, IHC, IF
NP Polyclonal Antibody	A55006	Influenza B virus	ELISA
NP Polyclonal Antibody	A55471	Influenza A virus	ELISA
NP Polyclonal Antibody	A56607	Influenza B virus	ELISA
NPBWR2 Polyclonal Antibody	A63038	Human	ELISA, IF
NPC2 Polyclonal Antibody	A56174	Dog	ELISA

Product Name	Cat. #	Species Reactivity	Applications
NPC2 Polyclonal Antibody	A60062	Human	ELISA, IHC, IF
NPHP1 Polyclonal Antibody	A53337	Human	ELISA, IHC, IF
NPHP4 Polyclonal Antibody	A68639	Human	ELISA, IHC, IF
NPHS1 Polyclonal Antibody	A56386	Human	ELISA, IHC
NPM2 Polyclonal Antibody	A65914	Human	ELISA, IHC
NPPB Polyclonal Antibody	A50644	Human, Mouse	ELISA, WB, IHC, IF
NPPB Polyclonal Antibody	A51533	Human	ELISA, WB, IHC
NPPB Polyclonal Antibody	A52796	Human	ELISA
NPPB Polyclonal Antibody	A54268	Human	ELISA
NPR1 Polyclonal Antibody	A56595	Human	ELISA
NPR2 Polyclonal Antibody	A56366	Human	ELISA
Npr3 Polyclonal Antibody	A63042	Rat	ELISA
NPY4R Polyclonal Antibody	A63210	Human	ELISA, IHC
Nqo2 Polyclonal Antibody	A50973	Rat, Mouse	ELISA, WB
NROB2 Polyclonal Antibody	A60066	Human	ELISA
NR1H3 Polyclonal Antibody	A55343	Human, Mouse	ELISA, WB, IHC
NR1H4 Polyclonal Antibody	A55491	Human, Mouse	ELISA, WB, IHC
NR1I3 Polyclonal Antibody	A55082	Human	ELISA, IHC
NR2C2AP Polyclonal Antibody	A63046	Human	ELISA, WB, IHC
NRARP Polyclonal Antibody	A63050	Human	ELISA, IHC
NRBP2 Polyclonal Antibody	A67610	Human	ELISA, IHC
NRCAM Polyclonal Antibody	A69635	Human	ELISA, IHC, IF
nrdA Polyclonal Antibody	A50813	Escherichia coli	ELISA
nrdB Polyclonal Antibody	A57401	Escherichia coli	ELISA
NRDC Polyclonal Antibody	A68643	Human	ELISA, WB, IF
nrdD Polyclonal Antibody	A53409	Escherichia coli	ELISA
NREP Polyclonal Antibody	A68871	Human	ELISA, IHC, IF
NRG2 Polyclonal Antibody	A60070	Human	ELISA, WB, IHC, IF
NRG4 Polyclonal Antibody	A64124	Human	ELISA, IHC, IF
NRIP2 Polyclonal Antibody	A67654	Human, Rat	ELISA, WB, IHC
NRIP3 Polyclonal Antibody	A60074	Human	ELISA, IHC
NRN1 Polyclonal Antibody	A55483	Human	ELISA, IHC
NRN1L Polyclonal Antibody	A64986	Human	ELISA, IHC, IF
NRSN1 Polyclonal Antibody	A66146	Human, Mouse	ELISA, WB, IF
NRSN2 Polyclonal Antibody	A67818	Human, Mouse	ELISA, WB, IF
NRXN3 Polyclonal Antibody	A67586	Human, Mouse	ELISA, WB
NSD2 Polyclonal Antibody	A69771	Human	ELISA, IHC, IF
NSDHL Polyclonal Antibody	A69871	Human, Rat	ELISA, WB, IHC, IF
Nsf Polyclonal Antibody	A52093	Rat	ELISA
NSMCE2 Polyclonal Antibody	A60078	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
NSUN2 Polyclonal Antibody	A68895	Human, Mouse, Rat	ELISA, WB, IF
NSUN3 Polyclonal Antibody	A67378	Human	ELISA
NSUN4 Polyclonal Antibody	A60082	Human	ELISA, WB, IHC
NSUN7 Polyclonal Antibody	A66398	Human	ELISA, IHC
NT5C1A Polyclonal Antibody	A57760	Human	ELISA, IHC, IF
NT5C2 Polyclonal Antibody	A68799	Human, Mouse	ELISA, WB, IF
NTF4 Polyclonal Antibody	A53377	Human	ELISA, IHC, IF
NTH1 Polyclonal Antibody	A52424	Saccharomyces cerevisiae	ELISA
NTMT1 Polyclonal Antibody	A56062	Human	ELISA
NTS Polyclonal Antibody	A56835	Human, Mouse	ELISA, WB
NUBPL Polyclonal Antibody	A60086	Human	ELISA, WB, IHC
Nuclease P1 Polyclonal Antibody	A64318	Penicillium citrinum	ELISA
Nucleoprotein Polyclonal Antibody	A55014	Rabies virus	ELISA
Nucleoprotein Polyclonal Antibody	A55507	Rabies virus	ELISA
Nucleoprotein Polyclonal Antibody	A57828	Influenza A virus	ELISA
NUDCD2 Polyclonal Antibody	A66662	Human	ELISA
NUDCD3 Polyclonal Antibody	A60090	Human	ELISA, WB
NUDT11 Polyclonal Antibody	A60094	Human	ELISA, WB, IHC
NUDT13 Polyclonal Antibody	A65818	Human	ELISA
NUDT14 Polyclonal Antibody	A63054	Human	ELISA
NUDT21 Polyclonal Antibody	A63058	Human	ELISA, WB, IHC, IF
NUDT5 Polyclonal Antibody	A70331	Human	ELISA, IHC, IF
NUDT9 Polyclonal Antibody	A60098	Human, Mouse	ELISA, WB
NUMBL Polyclonal Antibody	A70355	Human	ELISA, IHC, IF
NUP153 Polyclonal Antibody	A56539	Human	ELISA, IHC
NUP210 Polyclonal Antibody	A54950	Human	ELISA, IHC, IF
NUP35 Polyclonal Antibody	A66522	Human	ELISA, WB
NUP37 Polyclonal Antibody	A60102	Human	ELISA
NUP43 Polyclonal Antibody	A66774	Human	ELISA, WB, IHC
NUP58 Polyclonal Antibody	A60106	Human, Mouse	ELISA, WB, IHC
NUP85 Polyclonal Antibody	A67538	Human, Mouse	ELISA, WB, IHC, IF
NUPL2 Polyclonal Antibody	A63062	Human	ELISA, IHC
NUTM2F Polyclonal Antibody	A62558	Human	ELISA
NXF3 Polyclonal Antibody	A55946	Human	ELISA, IHC, IF
NXF5 Polyclonal Antibody	A63066	Human	ELISA, WB, IHC, IF
NXNL2 Polyclonal Antibody	A60110	Human	ELISA, IHC
NXPE1 Polyclonal Antibody	A60114	Human	ELISA, WB, IHC
NXPE3 Polyclonal Antibody	A51137	Human	ELISA, IHC
NXPH2 Polyclonal Antibody	A57429	Human	ELISA, IHC
OAZ3 Polyclonal Antibody	A63070	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
OBP2A Polyclonal Antibody	A67526	Human	ELISA, IHC
OCA2 Polyclonal Antibody	A68795	Human, Rat	ELISA, WB
OCC1 Polyclonal Antibody	A69127	Human	ELISA, IHC
OCIAD1 Polyclonal Antibody	A60118	Human, Mouse	ELISA, WB, IHC
ODC1 Polyclonal Antibody	A52669	Human	ELISA, WB, IHC, IF
ODF1 Polyclonal Antibody	A64654	Human	ELISA
ODF3L1 Polyclonal Antibody	A66122	Human	ELISA, WB, IHC
OFCC1 Polyclonal Antibody	A56659	Human, Mouse	ELISA, WB
OGDH Polyclonal Antibody	A55210	Human	ELISA, IHC
OGFOD2 Polyclonal Antibody	A65566	Human, Mouse	ELISA, WB
OGG1 Polyclonal Antibody	A60122	Human	ELISA, WB, IHC
OGN Polyclonal Antibody	A57045	Human, Rat	ELISA, WB, IHC
OGT Polyclonal Antibody	A56310	Human	ELISA, IHC, IF
OIH Polyclonal Antibody	A57716	Gallus gallus	ELISA
OIP5 Polyclonal Antibody	A63074	Human	ELISA
OLA1 Polyclonal Antibody	A54448	Human, Mouse	ELISA, WB, IHC, IF
OLFM4 Polyclonal Antibody	A68987	Human, Rat	ELISA, WB, IHC
OLFML1 Polyclonal Antibody	A65394	Human, Mouse	ELISA, WB, IHC, IF
OLFML2A Polyclonal Antibody	A65314	Human	ELISA, IHC
OLFML2B Polyclonal Antibody	A65334	Human, Mouse	ELISA, WB, IHC
OLR1 Polyclonal Antibody	A53032	Human	ELISA, IHC, IF
OMA1 Polyclonal Antibody	A67094	Human	ELISA, WB, IHC, IF
Omega-Gliadin Polyclonal Antibody	A57069	Triticum monococcum	ELISA
OMP Polyclonal Antibody	A69639	Human	ELISA, IF
ompA Polyclonal Antibody	A55519	Salmonella typhi	ELISA
ompA Polyclonal Antibody	A56338	Chlamydia trachomatis	ELISA
ompA Polyclonal Antibody	A56362	Chlamydophila pneumoniae	ELISA
ompA Polyclonal Antibody	A64052	Chlamydia pecorum	ELISA
ompC Polyclonal Antibody	A64454	Escherichia coli	ELISA
ONECUT1 Polyclonal Antibody	A69243	Human, Rat	ELISA, WB, IHC
OOSP2 Polyclonal Antibody	A65814	Human	ELISA, IHC
OPALIN Polyclonal Antibody	A66322	Human	ELISA, IF
OPN5 Polyclonal Antibody	A65386	Human	ELISA, IHC, IF
OR10W1 Polyclonal Antibody	A66226	Human	ELISA
OR1D2 Polyclonal Antibody	A63078	Human	ELISA
OR2H1 Polyclonal Antibody	A67554	Human, Mouse	ELISA, WB, IHC, IF
OR2L13 Polyclonal Antibody	A66350	Human, Mouse	ELISA, WB, IF
OR3A1 Polyclonal Antibody	A63082	Human	ELISA
OR5L1 Polyclonal Antibody	A66642	Human	ELISA, IHC
ORC4 Polyclonal Antibody	A54848	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
ORF1 Polyclonal Antibody	A57205	Hepatitis E virus genotype 1	ELISA
ORF2 Polyclonal Antibody	A57077	Hepatitis E virus genotype 1	ELISA
ORF3 Polyclonal Antibody	A56398	Hepatitis E virus genotype 1	ELISA
ORM2 Polyclonal Antibody	A52225	Human, Mouse	ELISA, WB
ORMDL1 Polyclonal Antibody	A60126	Human	ELISA, IHC, IF
ORMDL2 Polyclonal Antibody	A60130	Human	ELISA, IHC, IF
ORMDL3 Polyclonal Antibody	A70163	Human	ELISA, IHC, IF
OSBPL2 Polyclonal Antibody	A60134	Human, Mouse	ELISA, WB, IHC
OSBPL6 Polyclonal Antibody	A67806	Human	ELISA, IHC
OSER1 Polyclonal Antibody	A67326	Human, Mouse	ELISA, WB, IHC
OSGIN2 Polyclonal Antibody	A60138	Human	ELISA, WB, IHC
OSM Polyclonal Antibody	A52278	Human	ELISA
OSR2 Polyclonal Antibody	A60142	Human	ELISA, WB, IHC
OTC Polyclonal Antibody	A57433	Human	ELISA, IHC
OTUB1 Polyclonal Antibody	A53953	Human	ELISA, WB, IHC
OTUD6B Polyclonal Antibody	A60146	Human	ELISA, WB, IHC
OTUD7B Polyclonal Antibody	A69943	Human	ELISA, IHC, IF
OTX2 Polyclonal Antibody	A55299	Human, Mouse, Zebrafish	ELISA, WB, IHC
Outer Capsid Glycoprotein VP7 Polyclonal Antibody	A57501	Rotavirus A	ELISA
Outer Capsid Protein VP4 Polyclonal Antibody	A57461	Rotavirus A	ELISA
OV16 Polyclonal Antibody	A56623	Onchocerca volvulus	ELISA
OVOL1 Polyclonal Antibody	A63086	Human	ELISA, WB, IHC
Ovomucoid Polyclonal Antibody	A57565	Chicken	ELISA
Ovotransferrin, Allergen Gal d III, Conalbumin, Serum Transferrin Polyclonal Antibody	A57611	Chicken	ELISA
OXCT1 Polyclonal Antibody	A54776	Human	ELISA, WB, IHC
OXNAD1 Polyclonal Antibody	A51093	Human, Mouse	ELISA, WB
OXR1 Polyclonal Antibody	A60150	Human	ELISA, WB, IHC
OXSM Polyclonal Antibody	A67410	Human	ELISA, IHC
P2RY13 Polyclonal Antibody	A67166	Human	ELISA, WB, IF
p46 Polyclonal Antibody	A56443	Mycoplasma hyopneumoniae	ELISA
p46 Polyclonal Antibody	A56450	Mycoplasma hyopneumoniae	ELISA
p46 Polyclonal Antibody	A64250	Mycoplasma hyopneumoniae	ELISA
P4HA3 Polyclonal Antibody	A57317	Human	ELISA, IHC
P4HB Polyclonal Antibody	A52628	Human, Mouse	ELISA, WB, IHC
P4hb Polyclonal Antibody	A55066	Mouse	ELISA
P4HTM Polyclonal Antibody	A51337	Human	ELISA
PA Polyclonal Antibody	A68242	Influenza A virus	ELISA
PA2G4 Polyclonal Antibody	A60154	Human	ELISA, WB, IHC
PACRGL Polyclonal Antibody	A60158	Human	ELISA
PACS2 Polyclonal Antibody	A69643	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
PADI1 Polyclonal Antibody	A68050	Human	ELISA
PADI2 Polyclonal Antibody	A56390	Human, Mouse, Rat	ELISA, WB, IHC, IF
PADI2 Polyclonal Antibody	A56402	Human, Mouse	ELISA, WB
PADI3 Polyclonal Antibody	A68054	Human	ELISA, IHC
PAF1 Polyclonal Antibody	A66750	Human, Mouse	ELISA, WB, IHC
PAFAH1B2 Polyclonal Antibody	A60162	Human	ELISA, IHC, IF
PAFAH1B3 Polyclonal Antibody	A68867	Human	ELISA, IHC, IF
PAG2 Polyclonal Antibody	A69915	Bovine	ELISA
PAGE1 Polyclonal Antibody	A60166	Human	ELISA, IHC
PAGE2B Polyclonal Antibody	A65126	Human	ELISA, IHC
PAH Polyclonal Antibody	A63090	Human, Mouse	ELISA, WB, IHC, IF
PAIP2 Polyclonal Antibody	A67138	Human, Mouse	ELISA, WB, IHC, IF
PAK4 Polyclonal Antibody	A55818	Human	ELISA, IHC
pal Polyclonal Antibody	A64382	Legionella pneumophila	ELISA
PAM Polyclonal Antibody	A68647	Human	ELISA, IHC, IF
PAM16 Polyclonal Antibody	A53765	Human	ELISA, WB
PANK1 Polyclonal Antibody	A60170	Human	ELISA, IHC, IF
PANX2 Polyclonal Antibody	A70071	Human, Mouse, Rat	ELISA, WB, IHC, IF
Paox Polyclonal Antibody	A66010	Mouse	ELISA
PAPLN Polyclonal Antibody	A63094	Human	ELISA
PAPPA Polyclonal Antibody	A53093	Human	ELISA, IHC
PAQR5 Polyclonal Antibody	A67994	Human	ELISA, WB, IHC, IF
PAQR7 Polyclonal Antibody	A65762	Human	ELISA, IF
PARD3 Polyclonal Antibody	A51053	Human	ELISA, IHC
PARK2 Polyclonal Antibody	A57333	Human	ELISA, IHC, IF
PARK7 Polyclonal Antibody	A53957	Human, Mouse, Rat	ELISA, WB, IHC
PARK7 Polyclonal Antibody	A64064	Human, Mouse	ELISA, WB, IHC, IF
PARN Polyclonal Antibody	A53761	Human	ELISA, WB, IHC, IF
Parvalbumin Beta Polyclonal Antibody	A55154	Gadus morhua	ELISA
PASD1 Polyclonal Antibody	A60174	Human	ELISA, WB
PASK Polyclonal Antibody	A70035	Human	ELISA, IF
PAX2 Polyclonal Antibody	A53210	Human	ELISA
PAX3 Polyclonal Antibody	A50404	Human	ELISA, WB, IHC, IF
PAXIP1 Polyclonal Antibody	A65670	Human	ELISA, IHC
PBLD Polyclonal Antibody	A60178	Human, Mouse	ELISA, WB, IHC
PBXIP1 Polyclonal Antibody	A60182	Human	ELISA, WB
PCBD1 Polyclonal Antibody	A63102	Human	ELISA, IHC, IF
PCBP1 Polyclonal Antibody	A64658	Human	ELISA
PCBP4 Polyclonal Antibody	A60186	Human	ELISA
PCCB Polyclonal Antibody	A69647	Human, Mouse, Rat	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
PCDH11X Polyclonal Antibody	A70219	Human	ELISA, IHC, IF
PCDH20 Polyclonal Antibody	A66378	Human	ELISA
PCED1A Polyclonal Antibody	A67366	Human	ELISA, IHC
PCGF1 Polyclonal Antibody	A70195	Human	ELISA, IHC, IF
PCIF1 Polyclonal Antibody	A60190	Human	ELISA, IHC
PCK1 Polyclonal Antibody	A52968	Human	ELISA, IHC, IF
PCK2 Polyclonal Antibody	A64754	Human, Mouse	ELISA, WB
PCOLCE2 Polyclonal Antibody	A60194	Human	ELISA, WB, IHC, IF
PCP4 Polyclonal Antibody	A63106	Human	ELISA
PCSK2 Polyclonal Antibody	A68651	Human, Mouse	ELISA, WB, IHC
PCSK5 Polyclonal Antibody	A63110	Human	ELISA, IF
Pcsk5 Polyclonal Antibody	A63111	Mouse	ELISA
PCSK7 Polyclonal Antibody	A69651	Human	ELISA, IHC
PCTP Polyclonal Antibody	A63118	Human	ELISA, WB, IHC, IF
PDCD1 Polyclonal Antibody	A60198	Human, Mouse	ELISA, WB, IHC, IF
PDCD10 Polyclonal Antibody	A57153	Human	ELISA
PDCD1LG2 Polyclonal Antibody	A63122	Human	ELISA
PDCD6 Polyclonal Antibody	A51813	Human	ELISA, WB
PDCL2 Polyclonal Antibody	A60202	Human, Rat	ELISA, WB, IHC, IF
PDE6D Polyclonal Antibody	A60206	Human, Mouse	ELISA, WB, IHC
PDE6H Polyclonal Antibody	A64702	Human	ELISA
PDE9A Polyclonal Antibody	A68815	Human	ELISA, WB, IHC, IF
PDGFRB Polyclonal Antibody	A60210	Human	ELISA, IHC, IF
PDHA1 Polyclonal Antibody	A60214	Human	ELISA, IHC, IF
Pdia2 Polyclonal Antibody	A56186	Mouse, Human	ELISA, WB
PDIA3 Polyclonal Antibody	A51909	Human	ELISA
PDIA4 Polyclonal Antibody	A63126	Human, Mouse	ELISA, WB, IHC
PDIA5 Polyclonal Antibody	A60218	Human	ELISA, WB, IHC
PDIK1L Polyclonal Antibody	A60222	Human	ELISA, IHC
PDLIM4 Polyclonal Antibody	A68655	Human	ELISA, WB
PDP1 Polyclonal Antibody	A70279	Human	ELISA, WB, IHC
PDP2 Polyclonal Antibody	A55010	Human	ELISA
PDRG1 Polyclonal Antibody	A56939	Human	ELISA, IHC, IF
PDSS2 Polyclonal Antibody	A68999	Human	ELISA, IHC
PDXDC1 Polyclonal Antibody	A68979	Human	ELISA, IF
PDZD11 Polyclonal Antibody	A55371	Human	ELISA, IHC
PDZK1 Polyclonal Antibody	A54413	Human	ELISA
PECR Polyclonal Antibody	A60226	Human	ELISA
Pectate Lyase 1 Polyclonal Antibody	A50724	Juniperus ashei	ELISA
Pectate Lyase 1 Polyclonal Antibody	A51205	Chamaecyparis obtusa	ELISA



Product Name	Cat. #	Species Reactivity	Applications
Pectate Lyase 5 Polyclonal Antibody	A50788	Ambrosia artemisiifolia	ELISA
PEG10 Polyclonal Antibody	A54575	Human	ELISA, IHC
PELO Polyclonal Antibody	A60230	Human	ELISA, IHC
PEX11B Polyclonal Antibody	A60234	Human	ELISA, WB, IHC, IF
PEX13 Polyclonal Antibody	A69111	Human, Rat	ELISA, WB
PEX26 Polyclonal Antibody	A69959	Human	ELISA, WB, IHC
PF14_0076 Polyclonal Antibody	A65450	Plasmodium falciparum	ELISA
PFDN1 Polyclonal Antibody	A51489	Human	ELISA, IHC, IF
PFDN2 Polyclonal Antibody	A60238	Human	ELISA, WB, IHC, IF
PFDN5 Polyclonal Antibody	A51513	Human	ELISA, IHC, IF
PFDN5 Polyclonal Antibody	A51557	Human	ELISA, WB
PFKFB4 Polyclonal Antibody	A68831	Human	ELISA, IF
PFN1 Polyclonal Antibody	A51861	Human	ELISA, WB
PGA3 Polyclonal Antibody	A57701	Human	ELISA
PGC Polyclonal Antibody	A53492	Human, Mouse	ELISA, WB
PGD Polyclonal Antibody	A52692	Human	ELISA, WB
PGF Polyclonal Antibody	A60246	Human	ELISA
PGK1 Polyclonal Antibody	A51628	Human	ELISA, WB
PGK2 Polyclonal Antibody	A60250	Human	ELISA, WB
PGLYRP2 Polyclonal Antibody	A54408	Human	ELISA, IHC
PGLYRP4 Polyclonal Antibody	A63130	Human	ELISA, IHC
PGM1 Polyclonal Antibody	A50408	Human	ELISA, WB, IHC, IF
PGM2L1 Polyclonal Antibody	A65694	Human	ELISA, WB, IHC
PGR Polyclonal Antibody	A54696	Human	ELISA
PGS1 Polyclonal Antibody	A64890	Human, Mouse	ELISA, WB, IHC
PHB Polyclonal Antibody	A52978	Human, Mouse	ELISA, WB, IHC, IP
PHF10 Polyclonal Antibody	A60254	Human	ELISA, WB, IHC, IF
PHF13 Polyclonal Antibody	A63134	Human	ELISA, IHC
PHF5A Polyclonal Antibody	A53312	Human	ELISA, IHC, IF
PHF6 Polyclonal Antibody	A57345	Human	ELISA, WB, IHC
PHGDH Polyclonal Antibody	A51657	Human	ELISA, WB, IHC, IF
PHGDH Polyclonal Antibody	A53941	Human	ELISA, WB, IHC
Phosphoprotein Polyclonal Antibody	A54523	Rabies virus	ELISA, WB
PHOX2A Polyclonal Antibody	A68659	Human	ELISA, IHC
PHPT1 Polyclonal Antibody	A53656	Human	ELISA, WB, IHC, IF
phyA Polyclonal Antibody	A64346	Aspergillus niger	ELISA
PI3 Polyclonal Antibody	A60258	Human	ELISA, IHC, IF
PI-3 Polyclonal Antibody	A64282	Ascaris suum	ELISA
PI4KAP1 Polyclonal Antibody	A66206	Human, Mouse	ELISA, WB
PIAS2 Polyclonal Antibody	A50704	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
PICALM Polyclonal Antibody	A69903	Human	ELISA, IHC, IF
PICK1 Polyclonal Antibody	A55770	Human, Mouse	ELISA, WB, IHC
PID1 Polyclonal Antibody	A65942	Human	ELISA, IHC, IF
PIGC Polyclonal Antibody	A60262	Human	ELISA, IF
Pigeon Hemoglobin Polyclonal Antibody	A57600	Pigeon	ELISA
Pigeon Serum Albumin Polyclonal Antibody	A57528	Pigeon	ELISA
PIGH Polyclonal Antibody	A64706	Human, Mouse	ELISA, WB
PIGL Polyclonal Antibody	A68218	Human	ELISA, IHC
PIGO Polyclonal Antibody	A60266	Human	ELISA, IHC, IF
PIGR Polyclonal Antibody	A53372	Bovine	ELISA, WB
PIGS Polyclonal Antibody	A66826	Human	ELISA, IF
PIGV Polyclonal Antibody	A70227	Human	ELISA, WB, IHC, IF
PIGW Polyclonal Antibody	A69975	Human, Rat	ELISA, WB
PIGX Polyclonal Antibody	A56855	Human, Mouse	ELISA, WB, IHC
PIGZ Polyclonal Antibody	A65878	Human	ELISA, IF
PIH1D1 Polyclonal Antibody	A60270	Human	ELISA, WB, IHC
PIH1D3 Polyclonal Antibody	A67598	Human	ELISA, IHC, IF
PIK3AP1 Polyclonal Antibody	A68967	Human	ELISA, IF
PIK3IP1 Polyclonal Antibody	A70135	Human, Mouse	ELISA, WB, IF
PIK3R1 Polyclonal Antibody	A60274	Human	ELISA, IHC
PIK3R5 Polyclonal Antibody	A51105	Human	ELISA, IHC
PILRA Polyclonal Antibody	A68106	Human	ELISA, IHC
PILRB Polyclonal Antibody	A68026	Human	ELISA, IHC, IF
PIN1 Polyclonal Antibody	A60278	Human	ELISA, IHC
PIN4 Polyclonal Antibody	A70183	Human, Mouse, Rat	ELISA, WB, IHC, IF
PIP Polyclonal Antibody	A55387	Human	ELISA, IHC
PISD Polyclonal Antibody	A55870	Human	ELISA
PITPNC1 Polyclonal Antibody	A60282	Human	ELISA, IHC
PITPNM1 Polyclonal Antibody	A51465	Human	ELISA
Pka-C1 Polyclonal Antibody	A64270	Drosophila melanogaster	ELISA
PKD1L2 Polyclonal Antibody	A60286	Human	ELISA, IHC, IF
PKD2L2 Polyclonal Antibody	A67762	Human	ELISA, IHC, IF
Pklr Polyclonal Antibody	A50412	Rat	ELISA, WB
PKLR Polyclonal Antibody	A63138	Human	ELISA
PKM Polyclonal Antibody	A54959	Human	ELISA, IHC, IF
PKM Polyclonal Antibody	A63142	Human	ELISA
PKMYT1 Polyclonal Antibody	A54305	Human	ELISA, IHC
PLA1A Polyclonal Antibody	A69655	Human	ELISA, WB, IHC
PLA2G16 Polyclonal Antibody	A60290	Human	ELISA, IHC
PLA2G16 Polyclonal Antibody	A60291	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
PLA2G2A Polyclonal Antibody	A52245	Human	ELISA, IHC
Pla2g2a Polyclonal Antibody	A54844	Rat	ELISA
PLA2G2D Polyclonal Antibody	A55351	Human	ELISA, IHC
PLA2G5 Polyclonal Antibody	A54328	Mouse	ELISA
Pla2g5 Polyclonal Antibody	A54668	Rat	ELISA
PLA2G7 Polyclonal Antibody	A52996	Human, Mouse	ELISA, WB, IHC
PLA2G7 Polyclonal Antibody	A55383	Human, Mouse	ELISA, WB, IHC
PLA2R1 Polyclonal Antibody	A57425	Human	ELISA, IHC
PLAC1 Polyclonal Antibody	A67394	Human	ELISA
PLAC9 Polyclonal Antibody	A65098	Human	ELISA, IF
PLAT Polyclonal Antibody	A56467	Human, Mouse	ELISA, WB, IHC
PLAU Polyclonal Antibody	A52158	Human, Mouse	ELISA, WB, IHC
Plau Polyclonal Antibody	A53710	Rat, Human	ELISA, WB
PLAUR Polyclonal Antibody	A50636	Human, Mouse	ELISA, WB
Plbd2 Polyclonal Antibody	A54761	Mouse	ELISA, WB
PLCD4 Polyclonal Antibody	A51269	Human	ELISA, IHC
PLCL2 Polyclonal Antibody	A60298	Human	ELISA, IHC
PLCXD1 Polyclonal Antibody	A67978	Human	ELISA, WB, IHC
PLCZ1 Polyclonal Antibody	A65778	Human, Mouse	ELISA, WB
PLD Polyclonal Antibody	A56571	Loxosceles amazonica	ELISA
PLD4 Polyclonal Antibody	A51089	Human	ELISA, WB, IHC, IF
PLD5 Polyclonal Antibody	A55607	Human	ELISA, IHC
PLEC Polyclonal Antibody	A69883	Human	ELISA, IHC, IF
PLEKHA8 Polyclonal Antibody	A60302	Human, Mouse	ELISA, WB, IHC
PLEKHB2 Polyclonal Antibody	A56795	Human	ELISA, IHC
PLEKHJ1 Polyclonal Antibody	A67406	Human	ELISA, IHC
PLEKHS1 Polyclonal Antibody	A60306	Human	ELISA, WB, IHC
Plet1 Polyclonal Antibody	A60310	Mouse, Human	ELISA, WB, IHC
PLGRKT Polyclonal Antibody	A67582	Human, Mouse	ELISA, WB, IHC, IF
PLK1 Polyclonal Antibody	A63146	Human	ELISA, WB, IHC
PLLP Polyclonal Antibody	A68154	Human, Mouse	ELISA, WB, IF
PLN Polyclonal Antibody	A60314	Human	ELISA, IHC
PLOD2 Polyclonal Antibody	A63150	Human	ELISA, WB
pIp Polyclonal Antibody	A57301	Oncorhynchus mykiss	ELISA
PLP1 Polyclonal Antibody	A63154	Human	ELISA, WB, IHC, IF
PLPP2 Polyclonal Antibody	A60362	Human	ELISA, WB, IHC, IF
PLPP3 Polyclonal Antibody	A63190	Human, Mouse	ELISA, IHC, IF
PLTP Polyclonal Antibody	A53320	Human	ELISA, IHC
PLXNA1 Polyclonal Antibody	A52556	Human	ELISA, IHC
PLXNA4 Polyclonal Antibody	A67218	Human	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
ply Polyclonal Antibody	A50732	Streptococcus pneumoniae serotype 4	ELISA
PMP36 Polyclonal Antibody	A64394	Pig	ELISA
PMEL Polyclonal Antibody	A56587	Human	ELISA, WB
PMM1 Polyclonal Antibody	A50416	Human, Mouse	ELISA, WB, IHC
PMP2 Polyclonal Antibody	A51589	Human	ELISA, WB
PMP2 Polyclonal Antibody	A64262	Horse	ELISA
PMP22 Polyclonal Antibody	A60318	Human	ELISA, IF
PMRT15 Polyclonal Antibody	A69043	Arabidopsis thaliana	ELISA
PNLIPRP3 Polyclonal Antibody	A63158	Human, Mouse	ELISA, WB, IHC
PNMA1 Polyclonal Antibody	A60322	Human, Mouse	ELISA, WB, IHC
PNMA5 Polyclonal Antibody	A55331	Human	ELISA, IHC
PNO1 Polyclonal Antibody	A60326	Human	ELISA
PNP Polyclonal Antibody	A56551	Human	ELISA, WB
PNPLA1 Polyclonal Antibody	A66754	Human, Mouse	ELISA, WB, IHC, IF
PNPLA2 Polyclonal Antibody	A66446	Human, Mouse	ELISA, WB, IHC, IF
PNPLA5 Polyclonal Antibody	A69983	Human	ELISA, WB, IHC, IF
PNPLA8 Polyclonal Antibody	A69207	Human	ELISA, IHC, IF
PNPO Polyclonal Antibody	A69291	Human	ELISA, IHC, IF
PNRC2 Polyclonal Antibody	A60330	Human	ELISA, WB, IHC
pntB Polyclonal Antibody	A54459	Escherichia coli	ELISA
PODXL Polyclonal Antibody	A54939	Human	ELISA, WB, IHC, IF
POGK Polyclonal Antibody	A63162	Human	ELISA, IHC, IF
POGZ Polyclonal Antibody	A63166	Human	ELISA, IHC
poIA Polyclonal Antibody	A52038	Thermus aquaticus	ELISA, WB
POLA2 Polyclonal Antibody	A64674	Human	ELISA
POLB Polyclonal Antibody	A63170	Human	ELISA, WB, IHC
POLD2 Polyclonal Antibody	A69659	Human	ELISA, IHC
POLDIP2 Polyclonal Antibody	A68150	Human	ELISA
POLG Polyclonal Antibody	A69663	Human	ELISA, WB
POLG2 Polyclonal Antibody	A51361	Human	ELISA
POLK Polyclonal Antibody	A69295	Human, Mouse	ELISA, WB
POLR2C Polyclonal Antibody	A63174	Human, Mouse	ELISA, WB, IHC
POLR2D Polyclonal Antibody	A63178	Human	ELISA
POLR2E Polyclonal Antibody	A60334	Human	ELISA
POLR2I Polyclonal Antibody	A60338	Human	ELISA, IHC
POLR2J Polyclonal Antibody	A64402	Human	ELISA
POLR2K Polyclonal Antibody	A63182	Human	ELISA
POLR3A Polyclonal Antibody	A68316	Human	ELISA, IHC, CHIP
POLR3C Polyclonal Antibody	A55966	Human, Mouse	ELISA, WB
POLR3E Polyclonal Antibody	A60342	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
POLR3F Polyclonal Antibody	A60346	Human	ELISA, WB
POLR3K Polyclonal Antibody	A56102	Human	ELISA
Polygalacturonase Polyclonal Antibody	A50796	Cryptomeria japonica	ELISA
Polygalacturonase Polyclonal Antibody	A51005	Chamaecyparis obtusa	ELISA
POMC Polyclonal Antibody	A53537	Human	ELISA, IHC
POMC Polyclonal Antibody	A53538	Human	ELISA, WB
POMGNT1 Polyclonal Antibody	A60350	Human	ELISA, IHC
POMK Polyclonal Antibody	A67270	Human	ELISA, IHC
POMP Polyclonal Antibody	A63186	Human	ELISA
PON1 Polyclonal Antibody	A53284	Human, Mouse	ELISA, WB, IHC
PON3 Polyclonal Antibody	A68839	Human	ELISA, IHC
POP4 Polyclonal Antibody	A50420	Human	ELISA, IHC
POPDC3 Polyclonal Antibody	A69255	Human	ELISA, IF
POR Polyclonal Antibody	A57874	Human	ELISA, WB, IHC
Porcine Serum Albumin Polyclonal Antibody	A57557	Pig	ELISA
PORCN Polyclonal Antibody	A60354	Human, Mouse	ELISA, WB, IHC
pos5 Polyclonal Antibody	A68086	Schizosaccharomyces pombe	ELISA
POSTN Polyclonal Antibody	A54348	Human, Mouse	ELISA, WB, IHC, IF
POU2AF1 Polyclonal Antibody	A54276	Human	ELISA
POU5F1 Polyclonal Antibody	A55121	Human, Mouse	ELISA, WB, IHC
poxB Polyclonal Antibody	A52122	Escherichia coli	ELISA
PPAN Polyclonal Antibody	A60358	Human	ELISA
PPARG Polyclonal Antibody	A55407	Human	ELISA, IHC
PPBP Polyclonal Antibody	A52274	Human	ELISA
PPDK Polyclonal Antibody	A64358	Entamoeba histolytica	ELISA
PPEF1 Polyclonal Antibody	A60366	Human, Mouse	ELISA, WB, IHC
PPFIBP2 Polyclonal Antibody	A60370	Human	ELISA, WB
PPIA Polyclonal Antibody	A52820	Human, Mouse	ELISA, WB
PPIAL4G Polyclonal Antibody	A63194	Human, Mouse	ELISA, WB, IHC
PPIE Polyclonal Antibody	A70327	Human	ELISA, IHC
PPIH Polyclonal Antibody	A60374	Human	ELISA, WB, IHC
PPIL3 Polyclonal Antibody	A67254	Human	ELISA, WB, IHC
PPIL4 Polyclonal Antibody	A51169	Human, Mouse	ELISA, WB, IHC
PPM1K Polyclonal Antibody	A66354	Human, Mouse	ELISA, WB, IHC
PPM1M Polyclonal Antibody	A60378	Human	ELISA, IHC
PPME1 Polyclonal Antibody	A50424	Human, Mouse	ELISA, WB, IHC
PPOX Polyclonal Antibody	A63198	Human	ELISA, WB, IHC, IF, IP
PPP1CC Polyclonal Antibody	A63202	Human, Mouse	ELISA, WB, IHC
PPP1R11 Polyclonal Antibody	A51549	Human	ELISA, IHC
PPP1R14A Polyclonal Antibody	A66902	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
PPP1R16B Polyclonal Antibody	A66734	Human	ELISA, IHC
PPP1R3B Polyclonal Antibody	A65822	Human, Mouse, Rat	ELISA, WB, IHC
PPP1R7 Polyclonal Antibody	A60382	Human	ELISA, IHC
PPP2CA Polyclonal Antibody	A63206	Human, Mouse	ELISA, WB, IHC, IF
PPP2CB Polyclonal Antibody	A51581	Human	ELISA, WB, IHC, IF
PPP2R2B Polyclonal Antibody	A60386	Human	ELISA, WB
PPP2R3B Polyclonal Antibody	A68166	Human	ELISA
PPP2R4 Polyclonal Antibody	A60390	Human	ELISA, WB, IHC
PPP3R1 Polyclonal Antibody	A60394	Human	ELISA, WB
PPTC7 Polyclonal Antibody	A66530	Human	ELISA, IHC
PQLC2 Polyclonal Antibody	A60398	Human	ELISA, IF
PQLC3 Polyclonal Antibody	A60402	Human, Mouse	ELISA, WB, IHC, IF
PR10A Polyclonal Antibody	A64550	Coptis japonica	ELISA, WB
PRAC1 Polyclonal Antibody	A67110	Human	ELISA
PRAME Polyclonal Antibody	A56266	Human	ELISA, IHC
PRAMEF5 Polyclonal Antibody	A63214	Human	ELISA
PRCP Polyclonal Antibody	A68663	Human	ELISA, IF
PRDM14 Polyclonal Antibody	A67242	Human	ELISA, IHC
PRDM4 Polyclonal Antibody	A68034	Human, Mouse	ELISA, WB, IHC
PRDM5 Polyclonal Antibody	A70223	Human, Mouse	ELISA, WB
PRDX1 Polyclonal Antibody	A52589	Human	ELISA, WB, IHC, IF
PRDX4 Polyclonal Antibody	A60406	Human	ELISA, WB
Prealbumin Protein Polyclonal Antibody	A57604	Human	ELISA
PRELID2 Polyclonal Antibody	A66954	Human	ELISA
PRELID3A Polyclonal Antibody	A66722	Human	ELISA
PRELID3B Polyclonal Antibody	A61030	Human	ELISA, IHC
PRF Polyclonal Antibody	A57662	Toxoplasma gondii	ELISA
PRF1 Polyclonal Antibody	A52976	Human	ELISA
PRF1 Polyclonal Antibody	A63218	Mouse	ELISA
PRICKLE4 Polyclonal Antibody	A64882	Human	ELISA, IHC
PRIM1 Polyclonal Antibody	A63222	Human	ELISA, IF
PRKCH Polyclonal Antibody	A64176	Human, Mouse, Rat	ELISA, WB
PRKCSH Polyclonal Antibody	A51833	Human	ELISA, IHC
PRKRA Polyclonal Antibody	A68270	Human	ELISA, IHC
PRL Polyclonal Antibody	A60414	Human	ELISA
PRMT1 Polyclonal Antibody	A53937	Human, Mouse	ELISA, WB, IHC
PRMT6 Polyclonal Antibody	A60418	Human	ELISA, WB, IHC, IF
prn Polyclonal Antibody	A56499	Bordetella pertussis	ELISA
Prnp Polyclonal Antibody	A63226	Rat	ELISA
PROCR Polyclonal Antibody	A52584	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
PROK1 Polyclonal Antibody	A68667	Human	ELISA, IHC
Prok2 Polyclonal Antibody	A67650	Mouse	ELISA, WB
PROKR1 Polyclonal Antibody	A63230	Human	ELISA, IHC
PROS1 Polyclonal Antibody	A56286	Human	ELISA
PROSC Polyclonal Antibody	A60422	Human	ELISA, IHC
Protein Muc5ac Polyclonal Antibody	A57910	Mouse	ELISA
Protein P Polyclonal Antibody	A64298	Hepatitis B virus genotype A2 subtype adw2	ELISA
Protein X Polyclonal Antibody	A50877	Hepatitis B virus genotype D	ELISA
Protein X Polyclonal Antibody	A64286	Woodchuck hepatitis B virus	ELISA
Prothoracicotropic Hormone Polyclonal Antibody	A50772	Bombyx mori	ELISA
PRPF3 Polyclonal Antibody	A60426	Human	ELISA
PRPF8 Polyclonal Antibody	A69935	Human	ELISA, IF
PRPSAP1 Polyclonal Antibody	A60430	Human	ELISA, WB
PRR11 Polyclonal Antibody	A66458	Human, Mouse	ELISA, WB, IHC, IF
PRR3 Polyclonal Antibody	A56771	Human	ELISA
PRRC1 Polyclonal Antibody	A60434	Human	ELISA, IHC
PRRG3 Polyclonal Antibody	A67354	Human	ELISA, IF
Prss1 Polyclonal Antibody	A63234	Rat	ELISA
PRSS23 Polyclonal Antibody	A55918	Human	ELISA, IHC
Prss29 Polyclonal Antibody	A57509	Mouse	ELISA
PRSS3 Polyclonal Antibody	A52178	Human	ELISA, WB
PRSS35 Polyclonal Antibody	A66618	Human, Mouse	ELISA, WB, IHC
PRTFDC1 Polyclonal Antibody	A51377	Human	ELISA, IHC
PRY Polyclonal Antibody	A63238	Human	ELISA, IHC
PSAP Polyclonal Antibody	A53647	Human	ELISA, WB
Psat1 Polyclonal Antibody	A60438	Mouse	ELISA
PSAT1 Polyclonal Antibody	A68671	Human	ELISA, IHC, IF
PSD4 Polyclonal Antibody	A60442	Human	ELISA, IHC
PSG11 Polyclonal Antibody	A60446	Human	ELISA, WB
PSG3 Polyclonal Antibody	A60450	Human	ELISA, WB, IHC
PSG6 Polyclonal Antibody	A63242	Human	ELISA, WB
PSMA7 Polyclonal Antibody	A51517	Human, Mouse	ELISA, WB, IHC, IF
PSMB1 Polyclonal Antibody	A52597	Human	ELISA, WB
PSMB2 Polyclonal Antibody	A52680	Human, Mouse	ELISA, WB, IHC, IF
PSMB4 Polyclonal Antibody	A52715	Human	ELISA, WB, IHC, IF
PSMB7 Polyclonal Antibody	A52664	Human	ELISA, WB, IHC, IF
PSMC3 Polyclonal Antibody	A63246	Human	ELISA, IHC
PSMC6 Polyclonal Antibody	A60454	Human	ELISA
PSMD10 Polyclonal Antibody	A50428	Human	ELISA, IHC
PSMD11 Polyclonal Antibody	A51621	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
PSMD2 Polyclonal Antibody	A64758	Human	ELISA, WB, IHC, IF
PSMD7 Polyclonal Antibody	A60458	Human	ELISA, WB, IHC
PSME2 Polyclonal Antibody	A51485	Human	ELISA
PSME3 Polyclonal Antibody	A52301	Human	ELISA, IHC
PSMF1 Polyclonal Antibody	A57165	Human	ELISA, IHC
PSMG3 Polyclonal Antibody	A67174	Human	ELISA
PSTK Polyclonal Antibody	A60462	Human	ELISA, WB
PSTPIP2 Polyclonal Antibody	A60466	Human	ELISA, WB, IHC
pstS1 Polyclonal Antibody	A57705	Mycobacterium tuberculosis	ELISA
PTAFR Polyclonal Antibody	A63250	Human, Mouse	ELISA, WB, IHC, IF
PTCD2 Polyclonal Antibody	A66422	Human, Mouse	ELISA, WB, IHC
PTDSS1 Polyclonal Antibody	A51789	Human	ELISA, WB, IHC, IF
PTEN Polyclonal Antibody	A52436	Human	ELISA, IHC
PTGDS Polyclonal Antibody	A50432	Human	ELISA, IHC, IF
PTH Polyclonal Antibody	A52340	Human	ELISA, WB
PTH Polyclonal Antibody	A55810	Human	ELISA
pth Polyclonal Antibody	A64486	Escherichia coli	ELISA
PTH2R Polyclonal Antibody	A54886	Human	ELISA, IF
PTHLH Polyclonal Antibody	A54479	Human	ELISA, IHC
PTK2 Polyclonal Antibody	A52532	Chicken	ELISA
PTMA Polyclonal Antibody	A52464	Human	ELISA, IHC
PTP4A1 Polyclonal Antibody	A66810	Human	ELISA
PTP4A2 Polyclonal Antibody	A69843	Human	ELISA, IHC, IF
PTPMT1 Polyclonal Antibody	A55958	Human, Mouse, Rat	ELISA, WB, IHC
PTPN13 Polyclonal Antibody	A68835	Human	ELISA, IHC, IF
PTPN3 Polyclonal Antibody	A69667	Human	ELISA, IHC, IF
PTPN5 Polyclonal Antibody	A56098	Human	ELISA, WB, IHC
PTPN6 Polyclonal Antibody	A55619	Human	ELISA, WB, IHC, IF
PTPRN Polyclonal Antibody	A60470	Human	ELISA
PTPRR Polyclonal Antibody	A55271	Human	ELISA, IHC
PTPRZ1 Polyclonal Antibody	A56422	Human	ELISA, IHC, IF
PTPRZ1 Polyclonal Antibody	A56575	Human	ELISA, IHC
PTTG2 Polyclonal Antibody	A70275	Human	ELISA, IF
ptxA Polyclonal Antibody	A53718	Bordetella pertussis	ELISA, WB
PURA Polyclonal Antibody	A69671	Human	ELISA, WB
PVALB Polyclonal Antibody	A57009	Human, Mouse	ELISA, WB, IHC, IF
PVRIG Polyclonal Antibody	A65226	Human, Mouse	ELISA, WB, IF
PYC1 Polyclonal Antibody	A53675	Baker's yeast	ELISA
PYCR1 Polyclonal Antibody	A55723	Human	ELISA, IHC
PYCR2 Polyclonal Antibody	A60474	Human	ELISA, IHC



Product Name	Cat. #	Species Reactivity	Applications
PYGB Polyclonal Antibody	A51994	Human, Mouse	ELISA, WB, IHC
PYGL Polyclonal Antibody	A53440	Human	ELISA, WB, IHC
PYGO2 Polyclonal Antibody	A70307	Human	ELISA, IHC, IF
PYHIN1 Polyclonal Antibody	A65466	Human, Mouse	ELISA, WB
pyrF Polyclonal Antibody	A54629	Helicobacter pylori	ELISA
PYY Polyclonal Antibody	A54169	Human	ELISA, IHC
PZP Polyclonal Antibody	A53244	Human	ELISA
QDPR Polyclonal Antibody	A63254	Human, Mouse	ELISA, WB, IHC, IF
QPCT Polyclonal Antibody	A54495	Human, Mouse, Rat	ELISA, WB, IHC
QPCTL Polyclonal Antibody	A57341	Human	ELISA, IF
R107.2 Polyclonal Antibody	A64342	Caenorhabditis elegans	ELISA
R3HDML Polyclonal Antibody	A67490	Human	ELISA
RAB10 Polyclonal Antibody	A52568	Human	ELISA, IHC, IF
RAB11A Polyclonal Antibody	A56050	Human	ELISA, IHC
RAB11B Polyclonal Antibody	A60478	Human, Rat	ELISA, WB, IHC
RAB11FIP5 Polyclonal Antibody	A70199	Human	ELISA, IHC
RAB1A Polyclonal Antibody	A52624	Human	ELISA, WB
RAB20 Polyclonal Antibody	A67626	Human	ELISA, WB, IHC
RAB25 Polyclonal Antibody	A63258	Human, Mouse	ELISA, WB, IHC, IF
RAB27B Polyclonal Antibody	A50436	Human	ELISA
RAB2A Polyclonal Antibody	A55998	Human	ELISA
RAB30 Polyclonal Antibody	A50889	Human	ELISA, WB, IHC
RAB32 Polyclonal Antibody	A64730	Human, Mouse	ELISA, WB
RAB33A Polyclonal Antibody	A60482	Human	ELISA, IHC
RAB35 Polyclonal Antibody	A60486	Human	ELISA, IHC, IF
RAB37 Polyclonal Antibody	A60490	Human	ELISA, IHC
RAB38 Polyclonal Antibody	A60494	Human	ELISA
RAB39B Polyclonal Antibody	A66282	Human	ELISA, IHC, IF
RAB3B Polyclonal Antibody	A63262	Human, Mouse	ELISA, WB, IHC
RAB3C Polyclonal Antibody	A56779	Human, Mouse	ELISA, WB
RAB3IL1 Polyclonal Antibody	A60498	Human	ELISA
RAB40B Polyclonal Antibody	A60502	Human	ELISA, IHC
RAB4B Polyclonal Antibody	A63266	Human, Mouse	ELISA, WB
RAB5A Polyclonal Antibody	A56042	Human	ELISA, IHC
RAB6A Polyclonal Antibody	A57161	Human	ELISA, IHC
RAB6B Polyclonal Antibody	A60506	Human, Mouse	ELISA, WB, IHC
RAB6C Polyclonal Antibody	A67562	Human	ELISA, IHC
RAB8B Polyclonal Antibody	A66578	Human	ELISA, IHC
Rabbit Hemoglobin Polyclonal Antibody	A53073	Rabbit	ELISA
Rabbit Serum Albumin Polyclonal Antibody	A57592	Rabbit	ELISA

Product Name	Cat. #	Species Reactivity	Applications
RABEP2 Polyclonal Antibody	A67710	Human	ELISA, IHC
RABGEF1 Polyclonal Antibody	A51365	Human	ELISA, IHC
RABGGTA Polyclonal Antibody	A60510	Human, Mouse	ELISA, WB, IHC, IF
RABIF Polyclonal Antibody	A63270	Human, Mouse	ELISA, WB
RAC2 Polyclonal Antibody	A63274	Human	ELISA, IHC
RAD17 Polyclonal Antibody	A50440	Human, Mouse	ELISA, WB, IHC
RAD18 Polyclonal Antibody	A56282	Human	ELISA, WB, IHC
RAD23A Polyclonal Antibody	A53961	Human	ELISA, WB, IHC
RAD23B Polyclonal Antibody	A50444	Human	ELISA, WB, IHC
RAD54L2 Polyclonal Antibody	A69343	Human	ELISA, IF
RAD9B Polyclonal Antibody	A68947	Human, Mouse	ELISA, WB
RAF1 Polyclonal Antibody	A57169	Human	ELISA
RALGPS1 Polyclonal Antibody	A56382	Human	ELISA
RALY Polyclonal Antibody	A69315	Human, Rat	ELISA, WB, IHC, IF
RALYL Polyclonal Antibody	A60514	Human	ELISA, IHC
RANBP2 Polyclonal Antibody	A63278	Human	ELISA, IHC
RAP1A Polyclonal Antibody	A60518	Human	ELISA, IHC
RAP1B Polyclonal Antibody	A60522	Human, Mouse	ELISA, WB, IHC
RAPGEF5 Polyclonal Antibody	A60526	Human	ELISA, IHC
RARG Polyclonal Antibody	A68675	Human, Mouse	ELISA, WB, IHC
RARRES2 Polyclonal Antibody	A50664	Human	ELISA, IHC, IF
RARRES3 Polyclonal Antibody	A68114	Human	ELISA, IHC
RARS Polyclonal Antibody	A63282	Human	ELISA, WB, IHC, IF
RASA1 Polyclonal Antibody	A63286	Human, Mouse	ELISA, WB, IF
RASGEF1B Polyclonal Antibody	A64614	Human	ELISA, IHC
RASL10B Polyclonal Antibody	A66730	Human	ELISA, IHC
RASL12 Polyclonal Antibody	A67642	Human	ELISA
RASSF4 Polyclonal Antibody	A60530	Human	ELISA
RASSF7 Polyclonal Antibody	A70203	Human, Mouse	ELISA, WB, IF
Rat Serum Albumin Polyclonal Antibody	A57521	Rat	ELISA
RAVER2 Polyclonal Antibody	A55255	Human	ELISA, WB, IHC, IF
RBBP4 Polyclonal Antibody	A60534	Human, Mouse	ELISA, WB, IHC
RBBP7 Polyclonal Antibody	A64802	Human	ELISA, WB
RBBP9 Polyclonal Antibody	A60538	Human	ELISA, WB, IHC
RBCMT Polyclonal Antibody	A52010	Spinach	ELISA, WB
RBFOX2 Polyclonal Antibody	A68679	Human, Mouse	ELISA, WB, IHC, IF
RBFOX3 Polyclonal Antibody	A69835	Human	ELISA, IHC
RBKS Polyclonal Antibody	A60542	Human, Rat	ELISA, WB, IHC
RBL1 Polyclonal Antibody	A63290	Human, Mouse	ELISA, WB
RBM10 Polyclonal Antibody	A63294	Human, Mouse	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
RBM11 Polyclonal Antibody	A60546	Human	ELISA, IHC
RBM12 Polyclonal Antibody	A67310	Human	ELISA, IHC, IF
RBM22 Polyclonal Antibody	A67986	Human, Mouse	ELISA, WB, IHC
RBM3 Polyclonal Antibody	A50448	Human	ELISA, IHC
RBM41 Polyclonal Antibody	A60550	Human	ELISA, WB, IHC
RBM45 Polyclonal Antibody	A69047	Human, Rat	ELISA, WB, IF
RBM6 Polyclonal Antibody	A63298	Human	ELISA
RBM7 Polyclonal Antibody	A60554	Human	ELISA
RBMS2 Polyclonal Antibody	A60558	Human	ELISA, WB, IHC
RBMS3 Polyclonal Antibody	A68983	Human	ELISA, IF
RBMY1A1 Polyclonal Antibody	A57856	Human, Mouse, Rat	ELISA, WB, IHC, IF
RBMY1F Polyclonal Antibody	A50905	Human	ELISA, WB, IHC
RBP4 Polyclonal Antibody	A54045	Human	ELISA, WB, IHC
RBP4 Polyclonal Antibody	A54261	Human	ELISA
RBP4 Polyclonal Antibody	A57898	Human	ELISA, IHC
RBP4 Polyclonal Antibody	A64100	Human	ELISA
RBP5 Polyclonal Antibody	A60562	Human, Mouse	ELISA, WB
RBPMS Polyclonal Antibody	A60566	Human	ELISA, IHC, IF
RBPMS Polyclonal Antibody	A66274	Human	ELISA, IHC
rbsK Polyclonal Antibody	A52960	Escherichia coli	ELISA, WB
RBSN Polyclonal Antibody	A67822	Human	ELISA
rbtD Polyclonal Antibody	A52924	Enterobacter aerogenes	ELISA
RC3H2 Polyclonal Antibody	A67842	Human	ELISA
RCAN1 Polyclonal Antibody	A63302	Human	ELISA, IHC
RCBTB2 Polyclonal Antibody	A63306	Human, Rat	ELISA, WB, IHC, IF
RCOR3 Polyclonal Antibody	A63310	Human	ELISA, IHC, IF
RCVRN Polyclonal Antibody	A57281	Human, Mouse	ELISA, WB
RD3 Polyclonal Antibody	A65946	Human, Mouse	ELISA, WB, IHC
RDH11 Polyclonal Antibody	A55323	Human	ELISA, WB, IHC
RDH12 Polyclonal Antibody	A51117	Human	ELISA, IHC
recF Polyclonal Antibody	A64490	Escherichia coli	ELISA
recO Polyclonal Antibody	A64526	Escherichia coli	ELISA
recQ Polyclonal Antibody	A52364	Escherichia coli	ELISA, WB
recR Polyclonal Antibody	A64494	Escherichia coli	ELISA
REEP2 Polyclonal Antibody	A60570	Human	ELISA, IHC, IF
REEP3 Polyclonal Antibody	A65574	Human	ELISA, IHC, IF
REEP4 Polyclonal Antibody	A60574	Human	ELISA, WB, IHC, IF
REG1A Polyclonal Antibody	A63314	Human	ELISA, IHC
REG3A Polyclonal Antibody	A53824	Human	ELISA
Reg3a Polyclonal Antibody	A57505	Rat	ELISA

Product Name	Cat. #	Species Reactivity	Applications
Reg3b Polyclonal Antibody	A57245	Rat	ELISA
REG3G Polyclonal Antibody	A51001	Human	ELISA
Reg3g Polyclonal Antibody	A57073	Rat, Mouse	ELISA, WB
RELA Polyclonal Antibody	A68317	Human	ELISA, WB, IHC, ChIP
RELN Polyclonal Antibody	A57725	Human	ELISA, IHC
RELT Polyclonal Antibody	A52194	Human	ELISA, WB, IHC
Ren2 Polyclonal Antibody	A53129	Human, Mouse	ELISA, WB
repD Polyclonal Antibody	A64510	Staphylococcus aureus	ELISA
REPS1 Polyclonal Antibody	A70167	Human	ELISA, IHC, IF
RER1 Polyclonal Antibody	A60578	Human	ELISA, IHC, IF
RERG Polyclonal Antibody	A51153	Human, Mouse, Rat	ELISA, WB, IHC
RETN Polyclonal Antibody	A50632	Human	ELISA, IHC
REXO4 Polyclonal Antibody	A67238	Human	ELISA, IHC
RFESD Polyclonal Antibody	A66646	Human	ELISA, IHC
RFK Polyclonal Antibody	A66442	Human	ELISA, IHC, IF
RFPL3 Polyclonal Antibody	A60582	Human	ELISA, WB
RFPL4B Polyclonal Antibody	A63318	Human	ELISA, IHC, IF
RFTN2 Polyclonal Antibody	A65078	Human	ELISA, IHC, IF
RFX1 Polyclonal Antibody	A69675	Human	ELISA, WB, IHC
RFX2 Polyclonal Antibody	A68683	Human	ELISA, IF
RFX3 Polyclonal Antibody	A63322	Human	ELISA
RFX6 Polyclonal Antibody	A60586	Human	ELISA, IHC
RFXANK Polyclonal Antibody	A69679	Human	ELISA, WB, IHC, IF
RGMA Polyclonal Antibody	A66450	Human, Mouse	ELISA, WB, IHC
Rgn Polyclonal Antibody	A60590	Rat, Human	ELISA, WB
RGP1 Polyclonal Antibody	A60594	Human	ELISA
rgpA Polyclonal Antibody	A56206	Porphyromonas gingivalis	ELISA
RGS2 Polyclonal Antibody	A56997	Human, Mouse	ELISA, WB, IHC, IF
RGS22 Polyclonal Antibody	A63326	Human	ELISA
RGS4 Polyclonal Antibody	A63330	Human	ELISA, IHC, IF
RGS6 Polyclonal Antibody	A68687	Human, Rat	ELISA, WB, IHC, IF
RGS7 Polyclonal Antibody	A60598	Human	ELISA, IHC
RHBDL2 Polyclonal Antibody	A54699	Human, Mouse	ELISA, WB, IHC, IF
RHEBL1 Polyclonal Antibody	A60602	Human	ELISA, IHC
rho Polyclonal Antibody	A54197	Escherichia coli	ELISA
rho Polyclonal Antibody	A54875	Bacterium	ELISA
RHOA Polyclonal Antibody	A53873	Human	ELISA, IHC
RHOBTB3 Polyclonal Antibody	A63338	Human	ELISA, WB
RHOC Polyclonal Antibody	A60606	Human	ELISA, IHC
RHOD Polyclonal Antibody	A68691	Human	ELISA, IF

Product Name	Cat. #	Species Reactivity	Applications
RHOF Polyclonal Antibody	A67390	Human	ELISA, WB, IHC
RHOH Polyclonal Antibody	A64750	Human	ELISA, IHC, IF
RHOJ Polyclonal Antibody	A67494	Human	ELISA, IHC
RHOU Polyclonal Antibody	A68963	Human	ELISA, IHC, IF
RHOXF2 Polyclonal Antibody	A67170	Human	ELISA, IHC
RHPN1-AS1 Polyclonal Antibody	A67446	Human	ELISA, IHC
RIBC1 Polyclonal Antibody	A66486	Human	ELISA, IHC
RIBC2 Polyclonal Antibody	A67830	Human, Mouse	ELISA, WB, IHC
Riboflavin-Binding Protein Polyclonal Antibody	A55691	chicken	ELISA
Ribosome-Inactivating Protein Bryodin I Polyclonal Antibody	A64366	Bryonia dioica	ELISA
Ribosome-Inactivating Protein Charybdin Polyclonal Antibody	A64234	Drimia maritima	ELISA
Ribosome-Inactivating Protein Cucurmosin Polyclonal Antibody	A64222	Cucurbita moschata	ELISA, WB
Ribosome-inactivating Protein Gynostemmin Polyclonal Antibody	A65086	Gynostemma pentaphyllum	ELISA
Ribosome-Inactivating Protein Karasurin-C Polyclonal Antibody	A64306	Trichosanthes kirilowii	ELISA
Ribosome-Inactivating Protein Lychnin Polyclonal Antibody	A64226	Silene chalconica	ELISA
Ribosome-Inactivating Protein PD-L1/PD-L2 Polyclonal Antibody	A64216	Phytolacca dioica	ELISA
RIN1 Polyclonal Antibody	A68695	Human	ELISA, IF
RINT1 Polyclonal Antibody	A65474	Human	ELISA, IHC
RIPK2 Polyclonal Antibody	A54388	Human	ELISA, IHC
RIPK3 Polyclonal Antibody	A60610	Human	ELISA, IHC
RIT1 Polyclonal Antibody	A66806	Human	ELISA, IHC
RIT2 Polyclonal Antibody	A60614	Human	ELISA, IHC
RMDN1 Polyclonal Antibody	A66914	Human	ELISA, WB, IHC
RMDN3 Polyclonal Antibody	A56515	Human	ELISA, IHC
RMI1 Polyclonal Antibody	A67578	Human, Mouse	ELISA, WB
RMND5A Polyclonal Antibody	A55347	Human	ELISA, IHC
RMND5B Polyclonal Antibody	A60618	Human	ELISA, IHC
RNASE2 Polyclonal Antibody	A52888	Human	ELISA, IHC
RNASE3 Polyclonal Antibody	A64160	Human	ELISA
RNASE9 Polyclonal Antibody	A63342	Human	ELISA, IHC
RNASEL Polyclonal Antibody	A55090	Human	ELISA, IHC
RNF10 Polyclonal Antibody	A63346	Human, Mouse	ELISA, WB, IHC
RNF103 Polyclonal Antibody	A60622	Human	ELISA, IHC, IF
RNF128 Polyclonal Antibody	A66550	Human, Mouse	ELISA, WB, IF
RNF133 Polyclonal Antibody	A51193	Human	ELISA, WB, IHC, IF
RNF138 Polyclonal Antibody	A60626	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
RNF141 Polyclonal Antibody	A60630	Human	ELISA
RNF148 Polyclonal Antibody	A60634	Human	ELISA, WB, IHC, IF
RNF181 Polyclonal Antibody	A60638	Human	ELISA, IHC
RNF182 Polyclonal Antibody	A60642	Human	ELISA, IHC, IF
RNF183 Polyclonal Antibody	A60646	Human	ELISA, IHC, IF
RNF213 Polyclonal Antibody	A68923	Human	ELISA, IHC, IF
RNF24 Polyclonal Antibody	A60650	Human	ELISA, IHC, IF
RNF26 Polyclonal Antibody	A60654	Human	ELISA, WB, IHC, IF
RNF31 Polyclonal Antibody	A60658	Human	ELISA, IHC, IF
RNF39 Polyclonal Antibody	A68699	Human	ELISA, IF
RNF43 Polyclonal Antibody	A69683	Human	ELISA, IHC
RNF7 Polyclonal Antibody	A55050	Human	ELISA
rodA Polyclonal Antibody	A55802	Neosartorya fumigata	ELISA
ROGDI Polyclonal Antibody	A60666	Human	ELISA, WB, IHC
ROMO1 Polyclonal Antibody	A60670	Human	ELISA, IF
ROPN1 Polyclonal Antibody	A67962	Human	ELISA, IHC
ROPN1B Polyclonal Antibody	A54986	Human	ELISA, IHC
ROPN1L Polyclonal Antibody	A60674	Human	ELISA, IHC
ROR1 Polyclonal Antibody	A53536	Human	ELISA, IHC
RORC Polyclonal Antibody	A56258	Human	ELISA, IHC
RPA2 Polyclonal Antibody	A51461	Human	ELISA, WB, IHC, IF
RPA3 Polyclonal Antibody	A53417	Human	ELISA, WB
RPA3 Polyclonal Antibody	A57717	Human	ELISA, IHC
RPA4 Polyclonal Antibody	A54309	Human	ELISA
RPE Polyclonal Antibody	A66986	Human	ELISA, IHC, IF
RPE65 Polyclonal Antibody	A69911	Human, Rat	ELISA, WB, IF
RPGR Polyclonal Antibody	A60678	Human	ELISA, WB, IHC, IF
RPH3AL Polyclonal Antibody	A56523	Human, Rat	ELISA, WB
rpiA Polyclonal Antibody	A53512	Escherichia coli	ELISA, WB
RPL10L Polyclonal Antibody	A60682	Human, Mouse	ELISA, WB
RPL11 Polyclonal Antibody	A53841	Human, Mouse	ELISA, WB
RPL14 Polyclonal Antibody	A52992	Human	ELISA, IHC
RPL15 Polyclonal Antibody	A51741	Human	ELISA, WB
RPL17 Polyclonal Antibody	A51893	Human	ELISA, WB
RPL17 Polyclonal Antibody	A52704	Human	ELISA, WB
RPL26L1 Polyclonal Antibody	A68122	Human	ELISA, IHC
RPL27 Polyclonal Antibody	A52708	Human	ELISA, WB, IHC, IF
RPL28 Polyclonal Antibody	A53812	Human	ELISA, IHC, IF
RPL3 Polyclonal Antibody	A52836	Human	ELISA, IHC
RPL30 Polyclonal Antibody	A54978	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
RPL34 Polyclonal Antibody	A63350	Human	ELISA, IHC
RPL35A Polyclonal Antibody	A54938	Human	ELISA, IHC
RPL36AL Polyclonal Antibody	A51564	Human	ELISA, WB
RPL36AL Polyclonal Antibody	A51889	Human	ELISA, WB
RPL38 Polyclonal Antibody	A52728	Human	ELISA, WB, IF
RPL39L Polyclonal Antibody	A60686	Human	ELISA, IHC
RPL5 Polyclonal Antibody	A51577	Human	ELISA, WB, IHC
RPL7L1 Polyclonal Antibody	A65342	Human	ELISA, IHC
RPL8 Polyclonal Antibody	A52653	Human	ELISA, WB
RPLP0 Polyclonal Antibody	A57305	Human, Mouse, Rat	ELISA, WB, IHC, IF
RPLP1 Polyclonal Antibody	A52732	Human	ELISA, WB, IHC, IF
RPN1 Polyclonal Antibody	A54001	Human	ELISA, WB, IHC
RPN13 Polyclonal Antibody	A64586	Saccharomyces cerevisiae	ELISA
RPP14 Polyclonal Antibody	A60690	Human	ELISA, WB
RPP21 Polyclonal Antibody	A67954	Human	ELISA, WB
RPRD1A Polyclonal Antibody	A66930	Human	ELISA, WB
RPRD1B Polyclonal Antibody	A60694	Human	ELISA, WB, IHC
RPS10 Polyclonal Antibody	A52736	Human	ELISA, WB
RPS11 Polyclonal Antibody	A53308	Human	ELISA, WB, IHC
RPS12 Polyclonal Antibody	A52720	Human	ELISA, WB
RPS13 Polyclonal Antibody	A51576	Human	ELISA, WB, IHC
RPS14 Polyclonal Antibody	A51781	Human	ELISA
RPS16 Polyclonal Antibody	A52696	Human	ELISA, WB
RPS18 Polyclonal Antibody	A51757	Human	ELISA, WB
RPS19BP1 Polyclonal Antibody	A52812	Human	ELISA, WB, IHC
RPS24 Polyclonal Antibody	A60698	Human	ELISA, IHC
RPS25 Polyclonal Antibody	A51481	Human	ELISA
RPS25 Polyclonal Antibody	A51873	Human	ELISA
RPS26 Polyclonal Antibody	A63354	Human	ELISA, IHC
RPS28 Polyclonal Antibody	A60702	Human	ELISA, IHC
RPS29 Polyclonal Antibody	A69687	Human	ELISA, IF
RPS6KB2 Polyclonal Antibody	A53933	Human	ELISA, WB, IHC
RPS8 Polyclonal Antibody	A63358	Human, Mouse	ELISA, WB
RPS9 Polyclonal Antibody	A51877	Human	ELISA, WB
RPSA Polyclonal Antibody	A55158	Human, Mouse	ELISA, WB
rpsB Polyclonal Antibody	A51901	Escherichia coli	ELISA, WB
RPUSD4 Polyclonal Antibody	A60706	Human	ELISA, WB, IHC
RQCD1 Polyclonal Antibody	A68274	Human, Mouse	ELISA, WB, IHC
RRAGA Polyclonal Antibody	A51649	Human	ELISA, WB, IHC
RRM1 Polyclonal Antibody	A50452	Human	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
RRM1 Polyclonal Antibody	A60714	Human	ELISA, WB
RRM2 Polyclonal Antibody	A56054	Human	ELISA, WB
RS1 Polyclonal Antibody	A50456	Human	ELISA
RS1 Polyclonal Antibody	A57129	Human	ELISA, IHC
RSPH14 Polyclonal Antibody	A60718	Human	ELISA, WB, IHC
RSPH3 Polyclonal Antibody	A65802	Human	ELISA
RSPO3 Polyclonal Antibody	A60722	Human	ELISA, WB, IHC
RSPO3 Polyclonal Antibody	A67930	Human	ELISA, IHC
RSU1 Polyclonal Antibody	A50460	Human, Mouse	ELISA, WB
RTBDN Polyclonal Antibody	A60726	Human	ELISA
RTCA Polyclonal Antibody	A60730	Human	ELISA, IHC
rtcB Polyclonal Antibody	A55062	Escherichia coli	ELISA
RTKN2 Polyclonal Antibody	A60734	Human	ELISA, IHC
RTN4IP1 Polyclonal Antibody	A60738	Human	ELISA, IHC
RTN4RL1 Polyclonal Antibody	A65990	Human	ELISA, IHC, IF
RTP2 Polyclonal Antibody	A65270	Human, Mouse	ELISA, WB
RTP3 Polyclonal Antibody	A67902	Human, Mouse	ELISA, WB, IHC, IP
RTP4 Polyclonal Antibody	A60742	Human, Mouse	ELISA, WB, IHC
RUFY4 Polyclonal Antibody	A65538	Human	ELISA, IHC
RUNDC3A Polyclonal Antibody	A60746	Human	ELISA, WB, IHC
RUSC1 Polyclonal Antibody	A60750	Human	ELISA, IHC, IF
ruvA Polyclonal Antibody	A64498	Escherichia coli	ELISA
RUVBL1 Polyclonal Antibody	A68146	Human	ELISA, IHC
ruvC Polyclonal Antibody	A64530	Escherichia coli	ELISA
RWDD2A Polyclonal Antibody	A60754	Human, Mouse	ELISA, WB, IHC
RYR1 Polyclonal Antibody	A52748	Human	ELISA, IF
S100A1 Polyclonal Antibody	A55659	Human	ELISA, IHC
S100A10 Polyclonal Antibody	A63362	Human	ELISA, IHC, IF
S100A11 Polyclonal Antibody	A60758	Human	ELISA, IHC
S100A16 Polyclonal Antibody	A67002	Human	ELISA
S100A2 Polyclonal Antibody	A63366	Human	ELISA
S100A4 Polyclonal Antibody	A55109	Human	ELISA, WB, IHC
S100A6 Polyclonal Antibody	A55194	Human	ELISA, IHC
S100A8 Polyclonal Antibody	A53365	Human	ELISA
S100a8 Polyclonal Antibody	A56182	Mouse, Human	ELISA, IHC
S100a8 Polyclonal Antibody	A57117	Rat	ELISA
S100a8 Polyclonal Antibody	A57513	Mouse	ELISA
S100A9 Polyclonal Antibody	A53579	Human	ELISA, IHC
S100a9 Polyclonal Antibody	A55439	Mouse	ELISA
S100A9 Polyclonal Antibody	A60762	Human, Mouse	ELISA, WB



Product Name	Cat. #	Species Reactivity	Applications
S100B Polyclonal Antibody	A54129	Human	ELISA
S100g Polyclonal Antibody	A55219	Rat	ELISA
S100PBP Polyclonal Antibody	A60766	Human	ELISA, IHC
S1PR1 Polyclonal Antibody	A50464	Human, Mouse	ELISA, WB, IHC, IF
S1PR3 Polyclonal Antibody	A67154	Human	ELISA, WB, IF
SAA1 Polyclonal Antibody	A52372	Human	ELISA, WB, IHC
SAA1 Polyclonal Antibody	A55926	Dog	ELISA
sacIM Polyclonal Antibody	A57469	Streptomyces achromogenes	ELISA
SACM1L Polyclonal Antibody	A67402	Human	ELISA, IHC
SAG Polyclonal Antibody	A63370	Human	ELISA
SAMD12 Polyclonal Antibody	A66858	Human	ELISA, IHC
SAMD3 Polyclonal Antibody	A60770	Human	ELISA, IHC
SAMD7 Polyclonal Antibody	A65970	Human	ELISA, IHC
SAMHD1 Polyclonal Antibody	A53154	Mouse	ELISA, WB
SAMSN1 Polyclonal Antibody	A60774	Human	ELISA, IHC
SARAF Polyclonal Antibody	A66994	Human	ELISA
SARS Polyclonal Antibody	A53977	Human	ELISA, WB, IHC
SART3 Polyclonal Antibody	A60778	Human	ELISA, IHC
SAT1 Polyclonal Antibody	A53921	Human	ELISA, WB, IHC
SAT2 Polyclonal Antibody	A70043	Human	ELISA, IHC, IF
SBDS Polyclonal Antibody	A70351	Human	ELISA, IHC, IF
scaIM Polyclonal Antibody	A56851	Streptomyces caespitosus	ELISA
SCAMP1 Polyclonal Antibody	A68703	Human, Rat	ELISA, WB, IHC
SCAMP2 Polyclonal Antibody	A63374	Human	ELISA, IHC, IF
SCAMP3 Polyclonal Antibody	A54005	Human	ELISA, WB, IHC, IF
SCAMP4 Polyclonal Antibody	A60782	Human	ELISA, IHC, IF
SCAMP5 Polyclonal Antibody	A60786	Human	ELISA, IHC, IF
SCAND2P Polyclonal Antibody	A60790	Human	ELISA
SCAPER Polyclonal Antibody	A67230	Human	ELISA, IHC
SCARA3 Polyclonal Antibody	A65222	Human, Mouse	ELISA, WB, IHC
SCARB1 Polyclonal Antibody	A51125	Human	ELISA, IHC
SCFD2 Polyclonal Antibody	A60794	Human	ELISA, IHC
SCGB1C1 Polyclonal Antibody	A66534	Human	ELISA
SCGB2A1 Polyclonal Antibody	A63378	Human	ELISA
Scgb3a2 Polyclonal Antibody	A66802	Mouse	ELISA
SCGN Polyclonal Antibody	A63382	Human, Mouse	ELISA, WB, IHC, IF
SCLT1 Polyclonal Antibody	A63386	Human	ELISA, IHC
SCMH1 Polyclonal Antibody	A66818	Human	ELISA
scn Polyclonal Antibody	A67134	Staphylococcus aureus	ELISA
SCN10A Polyclonal Antibody	A68971	Human	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
SCN11A Polyclonal Antibody	A70303	Human	ELISA, IHC, IF
SCNM1 Polyclonal Antibody	A60798	Human	ELISA, IHC
SCNN1G Polyclonal Antibody	A60802	Human	ELISA, IHC, IF
SCOC Polyclonal Antibody	A63390	Human	ELISA, IHC
SCP Polyclonal Antibody	A57233	Epstein-Barr virus	ELISA
SCP2 Polyclonal Antibody	A54392	Human	ELISA, WB, IHC
SCRG1 Polyclonal Antibody	A60806	Human	ELISA, IHC
Scrg1 Polyclonal Antibody	A60807	Mouse	ELISA
SCRN3 Polyclonal Antibody	A60814	Human	ELISA
SCUBE2 Polyclonal Antibody	A69183	Human	ELISA, IHC
SCYL3 Polyclonal Antibody	A66050	Human	ELISA, IHC
Scytovirin Polyclonal Antibody	A64238	Scytonema varium	ELISA
SDC1 Polyclonal Antibody	A56799	Human	ELISA, IHC
SDCBP2 Polyclonal Antibody	A60818	Human	ELISA, IHC
SDCCAG3 Polyclonal Antibody	A69691	Human	ELISA, IHC, IF
SDF2 Polyclonal Antibody	A50468	Human	ELISA, IHC
SDF4 Polyclonal Antibody	A60822	Human	ELISA, WB
SDHA Polyclonal Antibody	A51569	Human	ELISA, WB, IHC, IF
SDHC Polyclonal Antibody	A60826	Human	ELISA, IHC, IF
SDHD Polyclonal Antibody	A54455	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC
SDSL Polyclonal Antibody	A63394	Human, Mouse	ELISA, WB, IHC
SEC11A Polyclonal Antibody	A60830	Human	ELISA, IHC
SEC11C Polyclonal Antibody	A57349	Human	ELISA, IHC
SEC14L3 Polyclonal Antibody	A67890	Human, Mouse	ELISA, WB, IHC
SEC14L4 Polyclonal Antibody	A67770	Human	ELISA, IHC
SEC16B Polyclonal Antibody	A67018	Human	ELISA, IHC
SEC22A Polyclonal Antibody	A60834	Human	ELISA, IHC, IF
SEC22B Polyclonal Antibody	A63398	Human	ELISA
SEC24C Polyclonal Antibody	A60838	Human	ELISA
SEC31A Polyclonal Antibody	A60842	Human	ELISA, IHC
SEC61G Polyclonal Antibody	A60846	Human	ELISA, IHC
SECTM1 Polyclonal Antibody	A60850	Human	ELISA, IHC, IF
SEL1L Polyclonal Antibody	A69187	Human	ELISA, IHC, IF
SELE Polyclonal Antibody	A63402	Sus scrofa	ELISA
SELM Polyclonal Antibody	A66566	Human	ELISA
Selt Polyclonal Antibody	A56326	Mouse	ELISA
SEMA5A Polyclonal Antibody	A64642	Human	ELISA, IHC
SEN8 Polyclonal Antibody	A56030	Human	ELISA
SEPP1 Polyclonal Antibody	A57909	Human, Mouse	ELISA, WB, IHC, IF
SEPSECS Polyclonal Antibody	A52925	Human	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
Septin 12 Polyclonal Antibody	A66046	Human, Mouse	ELISA, WB, IF
Septin 2 Polyclonal Antibody	A64718	Human	ELISA, IHC
SEPW1 Polyclonal Antibody	A63406	Human	ELISA, IHC
serB Polyclonal Antibody	A57225	Escherichia coli	ELISA
SERHL2 Polyclonal Antibody	A67190	Human	ELISA
Serine Protease Inhibitor Kazal-Type 2-Like Polyclonal Antibody	A56559	Cyprinus carpio	ELISA
Serotype 3 Fimbrial Subunit Polyclonal Antibody	A57848	Bordetella pertussis	ELISA
Serpina1 Polyclonal Antibody	A55133	Rat	ELISA
SERPINA1 Polyclonal Antibody	A64152	Human	ELISA
SERPINA11 Polyclonal Antibody	A69011	Human	ELISA, WB
Serpina3n Polyclonal Antibody	A66434	Mouse	ELISA
SERPINA6 Polyclonal Antibody	A54732	Human, Mouse, Rat	ELISA, WB, IHC
SERPINB1 Polyclonal Antibody	A55970	Human	ELISA, WB, IHC
SERPINB12 Polyclonal Antibody	A66318	Human	ELISA
SERPINB14 Polyclonal Antibody	A56374	Chicken	ELISA, WB
Serpib2 Polyclonal Antibody	A56222	Rat, Mouse	ELISA, WB
SERPINB3 Polyclonal Antibody	A55751	Human, Mouse	ELISA, WB, IHC
SERPINB4 Polyclonal Antibody	A68707	Human	ELISA, IHC, IF
SERPINB8 Polyclonal Antibody	A50472	Human	ELISA, WB, IHC
SERPINE1 Polyclonal Antibody	A52260	Human	ELISA, WB, IHC, IF
Serpinf1 Polyclonal Antibody	A50476	Rat	ELISA
Serpinh1 Polyclonal Antibody	A60854	Mouse, Human	ELISA, WB
SERTAD1 Polyclonal Antibody	A60858	Human	ELISA, WB, IHC
SES1 Polyclonal Antibody	A64458	Saccharomyces cerevisiae	ELISA
SETBP1 Polyclonal Antibody	A60862	Human	ELISA
SETD3 Polyclonal Antibody	A69963	Human	ELISA, IHC
SETD5 Polyclonal Antibody	A67470	Human	ELISA, IHC, IF
SETD9 Polyclonal Antibody	A66518	Human	ELISA, WB
SETX Polyclonal Antibody	A60870	Human	ELISA, IHC
SF3B6 Polyclonal Antibody	A68230	Human	ELISA, IHC
SFRP5 Polyclonal Antibody	A69695	Human	ELISA, IHC
SFT2D2 Polyclonal Antibody	A63410	Human, Mouse	ELISA, WB, IF
SFTA3 Polyclonal Antibody	A57057	Human	ELISA
SFTPB Polyclonal Antibody	A53476	Sus scrofa	ELISA
SFTPC Polyclonal Antibody	A55727	Bovine	ELISA
SFTPC Polyclonal Antibody	A60874	Human	ELISA, IHC
SFXN2 Polyclonal Antibody	A60878	Human	ELISA, IHC, IF
SFXN4 Polyclonal Antibody	A65434	Human	ELISA, WB, IHC, IF
SFXN5 Polyclonal Antibody	A66242	Human	ELISA
SGCA Polyclonal Antibody	A64830	Human	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
SGF29 Polyclonal Antibody	A58446	Human	ELISA, IHC
SGMS2 Polyclonal Antibody	A60882	Human	ELISA, IHC, IF
SGTB Polyclonal Antibody	A60886	Human	ELISA, IHC
SH3BGR Polyclonal Antibody	A60890	Human	ELISA
SH3BGR1 Polyclonal Antibody	A60894	Human	ELISA, IHC
SH3BGR13 Polyclonal Antibody	A67690	Human	ELISA, WB, IHC
SH3D19 Polyclonal Antibody	A65146	Human	ELISA, IHC
SH3GLB1 Polyclonal Antibody	A67590	Human	ELISA, IF
SHANK2 Polyclonal Antibody	A69327	Human	ELISA, IHC
SHARPIN Polyclonal Antibody	A69199	Human, Rat	ELISA, WB, IHC, IF
SHBG Polyclonal Antibody	A54447	Human	ELISA, WB
SHCBP1 Polyclonal Antibody	A66862	Human	ELISA
SHD Polyclonal Antibody	A66922	Human	ELISA, IHC
Sheep Serum Albumin Polyclonal Antibody	A57537	Sheep	ELISA
SHFM1 Polyclonal Antibody	A63414	Human	ELISA, IHC
SHISA5 Polyclonal Antibody	A60898	Human	ELISA, IHC
SHPRH Polyclonal Antibody	A69879	Human	ELISA, IHC, IF
SHROOM2 Polyclonal Antibody	A68887	Human	ELISA, IF
Sialic Acid-Binding Lectin Polyclonal Antibody	A56879	Rana japonica	ELISA
Sialidase-3 Polyclonal Antibody	A57815	Human	ELISA, WB
SIGLEC10 Polyclonal Antibody	A70095	Human	ELISA, IHC, IF
SIGLEC15 Polyclonal Antibody	A60902	Human, Mouse	ELISA, WB
SIGLEC6 Polyclonal Antibody	A60906	Human	ELISA, WB, IHC
SIL1 Polyclonal Antibody	A51333	Human, Mouse	ELISA, WB
SIM2 Polyclonal Antibody	A68879	Human	ELISA, IHC
SIMC1 Polyclonal Antibody	A60910	Human	ELISA
SIN3B Polyclonal Antibody	A68711	Human	ELISA, IHC, IF
SIPA1L2 Polyclonal Antibody	A60914	Human	ELISA, IHC
SIRPA Polyclonal Antibody	A50480	Human	ELISA, IHC
SIRPD Polyclonal Antibody	A56972	Human, Mouse	ELISA, WB
SIVA1 Polyclonal Antibody	A56891	Human	ELISA
SKA1 Polyclonal Antibody	A69107	Human	ELISA, IHC
SKA3 Polyclonal Antibody	A69051	Human	ELISA, IHC, IF
SKAP1 Polyclonal Antibody	A54280	Human	ELISA
SKP1 Polyclonal Antibody	A50484	Human, Mouse	ELISA, WB
SKP2 Polyclonal Antibody	A50885	Human	ELISA, WB, IHC
SLA Polyclonal Antibody	A60918	Human	ELISA, WB, IHC, IF
SLAIN1 Polyclonal Antibody	A69123	Human, Rat	ELISA, WB, IHC
SLAMF6 Polyclonal Antibody	A69699	Human, Mouse	ELISA, WB, IHC, IF
SLAMF8 Polyclonal Antibody	A63418	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
SLC10A2 Polyclonal Antibody	A64810	Human	ELISA, IHC, IF
SLC10A6 Polyclonal Antibody	A64950	Human	ELISA
SLC12A4 Polyclonal Antibody	A68126	Human	ELISA, IHC
SLC12A5 Polyclonal Antibody	A69179	Human	ELISA, IHC, IF
SLC12A9 Polyclonal Antibody	A60922	Human	ELISA, IF
SLC13A4 Polyclonal Antibody	A51397	Human	ELISA, WB, IHC, IF
SLC14A1 Polyclonal Antibody	A68847	Human	ELISA, IHC, IF
SLC17A3 Polyclonal Antibody	A63422	Human	ELISA, IHC, IF
SLC17A4 Polyclonal Antibody	A68202	Human	ELISA, IF
SLC19A2 Polyclonal Antibody	A69703	Human	ELISA, IHC
SLC1A6 Polyclonal Antibody	A63426	Human, Mouse	ELISA, WB, IF
SLC22A17 Polyclonal Antibody	A66418	Human	ELISA, IHC, IF
SLC22A18AS Polyclonal Antibody	A60926	Human	ELISA, IHC
SLC22A24 Polyclonal Antibody	A60930	Human	ELISA, IHC, IF
SLC22A9 Polyclonal Antibody	A66078	Human	ELISA, IHC, IF
SLC23A1 Polyclonal Antibody	A60934	Human, Mouse	ELISA, WB, IHC, IF
SLC23A3 Polyclonal Antibody	A60938	Human	ELISA, IHC, IF
SLC25A14 Polyclonal Antibody	A63430	Human	ELISA, IHC, IF
SLC25A18 Polyclonal Antibody	A67686	Human	ELISA, WB, IHC, IF
SLC25A25 Polyclonal Antibody	A60942	Human, Mouse	ELISA, WB, IHC, IF
SLC25A26 Polyclonal Antibody	A60946	Human	ELISA, IHC, IF
SLC25A32 Polyclonal Antibody	A67486	Human, Mouse	ELISA, WB
SLC25A33 Polyclonal Antibody	A60950	Human	ELISA, IHC, IF
SLC25A35 Polyclonal Antibody	A64970	Human	ELISA, IHC
SLC25A36 Polyclonal Antibody	A60954	Human	ELISA, WB, IHC, IF
SLC25A37 Polyclonal Antibody	A70251	Human	ELISA, IHC
SLC25A38 Polyclonal Antibody	A60958	Human	ELISA, IF
SLC25A4 Polyclonal Antibody	A60962	Human, Mouse	ELISA, WB, IHC
SLC25A40 Polyclonal Antibody	A60966	Human	ELISA, IHC, IF
SLC25A46 Polyclonal Antibody	A60970	Human, Mouse	ELISA, WB, IF
SLC25A51 Polyclonal Antibody	A60974	Human, Mouse	ELISA, WB, IF
SLC25A52 Polyclonal Antibody	A64942	Human	ELISA
SLC26A1 Polyclonal Antibody	A53925	Human	ELISA, IHC, IF
SLC26A4 Polyclonal Antibody	A69707	Human	ELISA, WB, IHC, IF
SLC26A6 Polyclonal Antibody	A69167	Human	ELISA, IHC, IF
SLC26A8 Polyclonal Antibody	A67030	Human	ELISA, IHC, IF
SLC27A5 Polyclonal Antibody	A54617	Human	ELISA
SLC28A1 Polyclonal Antibody	A63434	Human	ELISA, IHC, IF
SLC2A4 Polyclonal Antibody	A50488	Human, Mouse	ELISA, WB, IHC, IF
SLC2A4RG Polyclonal Antibody	A69711	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
SLC30A6 Polyclonal Antibody	A65430	Human	ELISA, WB, IHC, IF
SLC31A2 Polyclonal Antibody	A63438	Human	ELISA, IHC, IF
SLC34A3 Polyclonal Antibody	A70159	Human	ELISA, IHC, IF
SLC35B1 Polyclonal Antibody	A63442	Human	ELISA, WB, IF
SLC35B4 Polyclonal Antibody	A60978	Human	ELISA, IF
SLC35C1 Polyclonal Antibody	A60982	Human	ELISA
SLC35D2 Polyclonal Antibody	A65734	Human	ELISA, IHC
SLC35F2 Polyclonal Antibody	A66042	Human	ELISA
SLC35F5 Polyclonal Antibody	A60986	Human	ELISA, WB, IHC, IF
SLC35F6 Polyclonal Antibody	A60990	Human	ELISA, IHC, IF
SLC36A2 Polyclonal Antibody	A64990	Human	ELISA, IHC, IF
SLC36A3 Polyclonal Antibody	A64994	Human	ELISA, IHC, IF
SLC38A2 Polyclonal Antibody	A60994	Human	ELISA, WB, IHC, IF
SLC38A5 Polyclonal Antibody	A60998	Human	ELISA, IF
SLC38A6 Polyclonal Antibody	A66182	Human, Mouse	ELISA, WB, IF
SLC38A7 Polyclonal Antibody	A61002	Human	ELISA, IHC, IF
SLC39A11 Polyclonal Antibody	A61006	Human	ELISA, IF
Slc39a13 Polyclonal Antibody	A55643	Rat	ELISA
SLC41A2 Polyclonal Antibody	A53049	Human	ELISA, IHC, IF
SLC43A3 Polyclonal Antibody	A61010	Human	ELISA, IHC, IF
SLC44A4 Polyclonal Antibody	A50957	Human	ELISA, IF
SLC46A2 Polyclonal Antibody	A67454	Human, Mouse	ELISA, WB, IF
SLC46A3 Polyclonal Antibody	A65902	Human	ELISA, WB, IHC, IF
SLC4A8 Polyclonal Antibody	A61014	Human	ELISA, WB, IHC, IF
SLC52A1 Polyclonal Antibody	A67622	Human	ELISA, IHC, IF
SLC52A2 Polyclonal Antibody	A61018	Human	ELISA, IHC, IF
SLC5A10 Polyclonal Antibody	A63446	Human	ELISA, IF
SLC5A2 Polyclonal Antibody	A63450	Human	ELISA, IHC, IF
SLC5A8 Polyclonal Antibody	A70055	Human	ELISA, IF
SLC6A15 Polyclonal Antibody	A67566	Human	ELISA, IHC, IF
SLC6A16 Polyclonal Antibody	A61022	Human	ELISA, IHC, IF
SLC6A6 Polyclonal Antibody	A63454	Human	ELISA, WB, IHC, IF
SLC6A7 Polyclonal Antibody	A64072	Human	ELISA, IHC
SLC7A1 Polyclonal Antibody	A69715	Human	ELISA, IHC, IF
SLC7A14 Polyclonal Antibody	A61026	Human	ELISA, IHC
SLC7A2 Polyclonal Antibody	A68715	Human	ELISA, IHC
SLC7A3 Polyclonal Antibody	A67070	Human	ELISA, IF
SLC7A7 Polyclonal Antibody	A68118	Human, Mouse	ELISA, WB, IHC, IF
SLITRK3 Polyclonal Antibody	A63458	Human	ELISA
SLITRK4 Polyclonal Antibody	A66138	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
SLITRK6 Polyclonal Antibody	A67374	Human	ELISA, IHC, IF
SLPI Polyclonal Antibody	A51962	Human	ELISA, IHC, IF
SLURP1 Polyclonal Antibody	A63462	Human	ELISA, IHC, IF
SMAD1 Polyclonal Antibody	A61034	Human, Mouse	ELISA, WB, IHC
SMAD3 Polyclonal Antibody	A54649	Human	ELISA, WB
SMAD4 Polyclonal Antibody	A64766	Human	ELISA, WB, IHC, IP
SMAP1 Polyclonal Antibody	A61038	Human	ELISA, IHC
SMARCA2 Polyclonal Antibody	A68318	Human	ELISA, IHC, ChIP
SMARCB1 Polyclonal Antibody	A56154	Human	ELISA, IHC
SMC5 Polyclonal Antibody	A69999	Human	ELISA, WB
SMCO1 Polyclonal Antibody	A64842	Human	ELISA, IHC
SMCO3 Polyclonal Antibody	A62166	Human	ELISA, IF
SMCP Polyclonal Antibody	A53252	Human	ELISA, IHC
SMG6 Polyclonal Antibody	A65874	Human	ELISA, IHC, IF
SMIM14 Polyclonal Antibody	A61042	Human	ELISA, IHC, IF
SMIM19 Polyclonal Antibody	A61046	Human	ELISA, IHC, IF
SMNDC1 Polyclonal Antibody	A63466	Human	ELISA, IHC
Smox Polyclonal Antibody	A63470	Mouse	ELISA
Smox Polyclonal Antibody	A63474	Mouse, Human	ELISA, WB
SMPD2 Polyclonal Antibody	A68719	Human	ELISA, IHC, IF
SMPD3 Polyclonal Antibody	A67638	Human	ELISA, IHC
SMPDL3A Polyclonal Antibody	A61050	Human	ELISA, IHC
SMT3 Polyclonal Antibody	A50909	Saccharomyces cerevisiae	ELISA
SMT3 Polyclonal Antibody	A64698	Saccharomyces cerevisiae	ELISA
SMUG1 Polyclonal Antibody	A61054	Human	ELISA
SMURF2 Polyclonal Antibody	A61058	Human	ELISA, IHC
SNAI1 Polyclonal Antibody	A61062	Human	ELISA, IHC
SNAPC1 Polyclonal Antibody	A61066	Human	ELISA
SNCA Polyclonal Antibody	A54037	Human	ELISA, WB, IHC
Snca Polyclonal Antibody	A55267	Rat	ELISA
SNCA Polyclonal Antibody	A56511	Human	ELISA, IHC
SNCG Polyclonal Antibody	A56479	Human, Mouse	ELISA, WB, IHC
SNPH Polyclonal Antibody	A61070	Human	ELISA
SNRNP25 Polyclonal Antibody	A61074	Human	ELISA
Snrnp35 Polyclonal Antibody	A67810	Mouse, Human	ELISA, WB
SNRNP40 Polyclonal Antibody	A66286	Human	ELISA
SNRNP48 Polyclonal Antibody	A65502	Human	ELISA
SNRPA Polyclonal Antibody	A52652	Human, Mouse	ELISA, WB, IHC, IF
SNRPB2 Polyclonal Antibody	A50837	Human	ELISA, IHC
SNRPD1 Polyclonal Antibody	A63478	Human	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
SNRPD2 Polyclonal Antibody	A51653	Human	ELISA, WB
SNRPG Polyclonal Antibody	A53897	Human	ELISA, WB, IHC, IF
SNX12 Polyclonal Antibody	A70335	Human	ELISA, IHC, IF
SNX16 Polyclonal Antibody	A61078	Human	ELISA, IHC
SNX20 Polyclonal Antibody	A53539	Human, Mouse	ELISA, WB, IHC
SNX22 Polyclonal Antibody	A61082	Human	ELISA, IHC
SNX24 Polyclonal Antibody	A61086	Human	ELISA, IHC
SNX3 Polyclonal Antibody	A51509	Human	ELISA
SNX3 Polyclonal Antibody	A51881	Human	ELISA
SNX6 Polyclonal Antibody	A61090	Human	ELISA, WB, IHC
SOCS1 Polyclonal Antibody	A54436	Human	ELISA, IHC
Socs3 Polyclonal Antibody	A61094	Rat, Human	ELISA, WB
SOCS3 Polyclonal Antibody	A63482	Human	ELISA, IHC, IF
SOD1 Polyclonal Antibody	A52404	Human, Mouse	ELISA, WB, IF
Sod1 Polyclonal Antibody	A53702	Mouse, Human	ELISA, IHC
Sod1 Polyclonal Antibody	A54676	Mouse	ELISA
SOD2 Polyclonal Antibody	A50492	Human, Mouse	ELISA, WB, IHC
SOD3 Polyclonal Antibody	A61098	Human	ELISA
sodB Polyclonal Antibody	A66982	Neosartorya fumigata	ELISA
SOHLH2 Polyclonal Antibody	A61102	Human	ELISA, IHC
SORBS3 Polyclonal Antibody	A61106	Human	ELISA, WB, IHC
Sort1 Polyclonal Antibody	A56294	Rat	ELISA
SOSTDC1 Polyclonal Antibody	A53722	Human	ELISA
SOX2 Polyclonal Antibody	A54440	Human, Mouse	ELISA, WB, IHC
SOX21 Polyclonal Antibody	A69339	Human	ELISA, IF
SOX3 Polyclonal Antibody	A68723	Human	ELISA, WB, IHC
SOX8 Polyclonal Antibody	A69719	Human	ELISA, WB
SP6 Polyclonal Antibody	A64974	Human	ELISA, IHC
SPA17 Polyclonal Antibody	A52897	Human	ELISA
SPACA1 Polyclonal Antibody	A67290	Human	ELISA, IHC
SPACA9 Polyclonal Antibody	A66586	Human	ELISA
SPAG8 Polyclonal Antibody	A69155	Human	ELISA, IF
SPAM1 Polyclonal Antibody	A50496	Human	ELISA
SPANXC Polyclonal Antibody	A61110	Human	ELISA, IHC
SPATA19 Polyclonal Antibody	A65838	Human, Mouse	ELISA, WB
SPATA8 Polyclonal Antibody	A65710	Human	ELISA
SPATC1L Polyclonal Antibody	A67362	Human	ELISA, IHC, IF
SPATS1 Polyclonal Antibody	A65014	Human	ELISA, IHC
SPATS2L Polyclonal Antibody	A55974	Human, Mouse	ELISA, WB, IHC
SPATS2L Polyclonal Antibody	A61114	Human	ELISA, WB



Product Name	Cat. #	Species Reactivity	Applications
SPC25 Polyclonal Antibody	A67722	Human	ELISA, IHC
SPCS1 Polyclonal Antibody	A61118	Human	ELISA, IF
SPCS2 Polyclonal Antibody	A64714	Human	ELISA, WB, IHC, IF
SPCS3 Polyclonal Antibody	A63486	Human, Mouse	ELISA, WB, IHC
SPDL1 Polyclonal Antibody	A66698	Human	ELISA, WB, IHC
speA Polyclonal Antibody	A52213	Escherichia coli	ELISA
SPEM1 Polyclonal Antibody	A66366	Human	ELISA, IHC
SPERT Polyclonal Antibody	A61122	Human, Mouse, Rat	ELISA, WB
SPG11 Polyclonal Antibody	A66706	Human	ELISA, IF
SPG21 Polyclonal Antibody	A68002	Human	ELISA, IHC, IF
SPI1 Polyclonal Antibody	A54236	Human	ELISA
SPIC Polyclonal Antibody	A61126	Human, Mouse	ELISA, WB, IHC
Spike Glycoprotein Polyclonal Antibody	A56214	Human coronavirus OC43	ELISA
Spike Glycoprotein Polyclonal Antibody	A57201	Bovine coronavirus	ELISA
SPIN2A Polyclonal Antibody	A67146	Human	ELISA
SPIN3 Polyclonal Antibody	A65134	Human	ELISA, WB
SPIN4 Polyclonal Antibody	A65110	Human	ELISA, IHC, IF
SPINK6 Polyclonal Antibody	A68955	Human	ELISA, IHC
SPIRE2 Polyclonal Antibody	A51201	Human	ELISA, IHC, IF
SPOP Polyclonal Antibody	A61130	Human	ELISA, IHC
SPOPL Polyclonal Antibody	A69723	Human	ELISA, WB
SPP1 Polyclonal Antibody	A54033	Human	ELISA, IHC
Spp1 Polyclonal Antibody	A54808	Mouse	ELISA, WB
SPPL2A Polyclonal Antibody	A67058	Human	ELISA, WB, IHC, IF
SPRR1B Polyclonal Antibody	A63490	Human	ELISA, IHC
SPRR2A Polyclonal Antibody	A66690	Human	ELISA
SPRTN Polyclonal Antibody	A61134	Human	ELISA, IHC
SPRY1 Polyclonal Antibody	A61138	Human, Mouse	ELISA, WB, IHC
SPRY4 Polyclonal Antibody	A61142	Human	ELISA
SPT16 Polyclonal Antibody	A50512	Arabidopsis thaliana	ELISA
SPTBN1 Polyclonal Antibody	A68943	Human	ELISA, IF
SPTLC2 Polyclonal Antibody	A50500	Human	ELISA, WB, IHC, IF
SPTLC3 Polyclonal Antibody	A67314	Human, Mouse	ELISA, WB, IF
SPZ1 Polyclonal Antibody	A67926	Human, Mouse	ELISA, WB, IHC
SRFBP1 Polyclonal Antibody	A61146	Human	ELISA
SRGAP3 Polyclonal Antibody	A69727	Human	ELISA, IF
SRM Polyclonal Antibody	A53881	Human	ELISA, WB, IHC
SRP19 Polyclonal Antibody	A54760	Human	ELISA, WB, IHC
SRP19 Polyclonal Antibody	A63494	Human	ELISA, IHC, IF
SRP54 Polyclonal Antibody	A63498	Human, Mouse	ELISA, WB, IF

Product Name	Cat. #	Species Reactivity	Applications
SRPK3 Polyclonal Antibody	A68142	Human	ELISA, IHC
SRPRB Polyclonal Antibody	A51421	Human	ELISA, WB, IF
SRSF10 Polyclonal Antibody	A50873	Human	ELISA, IHC
SRSF9 Polyclonal Antibody	A51472	Human, Mouse	ELISA, WB
Sry Polyclonal Antibody	A53594	Mouse	ELISA
ssaA1 Polyclonal Antibody	A65046	Staphylococcus aureus	ELISA
SSB Polyclonal Antibody	A54756	Human, Mouse	ELISA, WB
ssb Polyclonal Antibody	A56727	Bacillus anthracis	ELISA
SSB Protein gp2.5 Polyclonal Antibody	A56230	Enterobacteria phage T7	ELISA
SSFA2 Polyclonal Antibody	A50504	Human	ELISA, IHC
SSH3 Polyclonal Antibody	A69075	Human, Rat	ELISA, WB, IHC, IF
SSMEM1 Polyclonal Antibody	A61150	Human	ELISA, IHC, IF
sspB Polyclonal Antibody	A54567	Staphylococcus aureus	ELISA
SSR3 Polyclonal Antibody	A61154	Human	ELISA, WB, IHC, IF
SSSCA1 Polyclonal Antibody	A56755	Human	ELISA, WB, IHC
SST Polyclonal Antibody	A63502	Human	ELISA, IHC
SST Polyclonal Antibody	A63503	Human	ELISA, IHC
SSU72 Polyclonal Antibody	A61158	Human	ELISA
SSX2 Polyclonal Antibody	A50893	Human	ELISA
ST3GAL2 Polyclonal Antibody	A50913	Human	ELISA
ST3GAL3 Polyclonal Antibody	A54288	Human, Mouse	ELISA, WB
ST3GAL4 Polyclonal Antibody	A69839	Human	ELISA, IHC
ST6GALNAC2 Polyclonal Antibody	A67786	Human	ELISA, WB
ST6GALNAC3 Polyclonal Antibody	A66514	Human	ELISA
ST6GALNAC6 Polyclonal Antibody	A61162	Human	ELISA, IHC
ST8SIA4 Polyclonal Antibody	A66266	Human	ELISA
ST8SIA5 Polyclonal Antibody	A63504	Human	ELISA, WB, IHC
STAC3 Polyclonal Antibody	A61166	Human, Rat	ELISA, WB
STAMBIP Polyclonal Antibody	A55623	Human	ELISA, WB, IHC
STAP1 Polyclonal Antibody	A61170	Human	ELISA
STAR Polyclonal Antibody	A61174	Human	ELISA, IHC
STARD10 Polyclonal Antibody	A51413	Human, Mouse	ELISA, WB
STARD7 Polyclonal Antibody	A53674	Human	ELISA, WB
STAT1 Polyclonal Antibody	A63508	Human	ELISA, IHC
STC1 Polyclonal Antibody	A61178	Human	ELISA, IHC
STEAP1 Polyclonal Antibody	A54181	Human	ELISA, WB, IHC, IF
STEAP3 Polyclonal Antibody	A61182	Human	ELISA, WB, IHC, IF
STIP1 Polyclonal Antibody	A54796	Human	ELISA, WB, IHC
STK17B Polyclonal Antibody	A63512	Human, Mouse	ELISA, WB, IHC
STK24 Polyclonal Antibody	A54344	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
STK25 Polyclonal Antibody	A54396	Human	ELISA, IHC
STK33 Polyclonal Antibody	A61186	Human, Mouse	ELISA, WB, IHC, IF
STK38L Polyclonal Antibody	A61190	Human	ELISA, WB, IHC
STOML1 Polyclonal Antibody	A67886	Human	ELISA, IF
STOX1 Polyclonal Antibody	A65730	Human	ELISA, WB, IHC
STPG2 Polyclonal Antibody	A66850	Human	ELISA, WB, IHC
STRA13 Polyclonal Antibody	A63516	Human	ELISA, IHC
STRA6 Polyclonal Antibody	A69191	Human	ELISA, IHC, IF
STRADB Polyclonal Antibody	A51325	Human	ELISA, WB, IHC
STRBP Polyclonal Antibody	A66830	Human	ELISA, WB, IF
Streptavidin Polyclonal Antibody	A50712	Streptomyces avidinii	ELISA
STT3A Polyclonal Antibody	A63520	Human	ELISA, IHC
STX10 Polyclonal Antibody	A61194	Human	ELISA, IHC
STX10 Polyclonal Antibody	A61195	Human	ELISA, IHC
STX19 Polyclonal Antibody	A66362	Human	ELISA, WB
STX4 Polyclonal Antibody	A61202	Human	ELISA
STX6 Polyclonal Antibody	A57693	Chicken, Mouse	ELISA, WB
STX6 Polyclonal Antibody	A57697	Human	ELISA, IHC
STX8 Polyclonal Antibody	A68098	Human	ELISA, IHC
stxB Polyclonal Antibody	A50720	Enterobacteria phage H19B	ELISA
stxB2 Polyclonal Antibody	A50821	Enterobacteria phage 933W	ELISA
STXBP6 Polyclonal Antibody	A51165	Human, Mouse	ELISA, WB, IHC, IF
SUB1 Polyclonal Antibody	A55129	Human	ELISA, IHC
SUCLG2 Polyclonal Antibody	A70099	Human	ELISA, WB, IHC
SUCNR1 Polyclonal Antibody	A63524	Human, Mouse	ELISA, WB, IF
sulA Polyclonal Antibody	A64534	Escherichia coli	ELISA
SULT1B1 Polyclonal Antibody	A50508	Human, Mouse	ELISA, WB, IHC
SUOX Polyclonal Antibody	A54789	Human	ELISA, WB, IHC
SUPT3H Polyclonal Antibody	A54487	Human, Mouse	ELISA, WB, IHC
surA Polyclonal Antibody	A57221	Escherichia coli	ELISA, WB
SUS1 Polyclonal Antibody	A64354	Saccharum officinarum	ELISA
SUSD4 Polyclonal Antibody	A61206	Human	ELISA, IHC
SUSD6 Polyclonal Antibody	A66438	Human	ELISA, IHC, IF
SV2B Polyclonal Antibody	A61210	Human, Mouse	ELISA, WB, IF
SVBP Polyclonal Antibody	A66482	Human	ELISA
SVOPL Polyclonal Antibody	A51041	Human, Mouse, Rat	ELISA, WB, IF
SVSP Polyclonal Antibody	A57814	Crotalus adamanteus	ELISA
SWAP70 Polyclonal Antibody	A61214	Human, Mouse	ELISA, WB
SYCN Polyclonal Antibody	A63528	Human	ELISA, IHC
SYF2 Polyclonal Antibody	A61218	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
SYN1 Polyclonal Antibody	A54720	Human, Mouse	ELISA, WB, IHC
SYNC Polyclonal Antibody	A67194	Human	ELISA
SYNCRIP Polyclonal Antibody	A51796	Human	ELISA, WB, IHC
SYNE2 Polyclonal Antibody	A70087	Human	ELISA, IHC, IF
SYNE4 Polyclonal Antibody	A66846	Human	ELISA
SYNJ1 Polyclonal Antibody	A69731	Human	ELISA, IHC, IF
SYNJ2BP Polyclonal Antibody	A63532	Human, Rat	ELISA, WB, IHC
SYNPR Polyclonal Antibody	A61222	Human, Mouse, Rat	ELISA, WB, IHC, IF
SYT1 Polyclonal Antibody	A52536	Human	ELISA
SYT10 Polyclonal Antibody	A65598	Human	ELISA, IHC
SYTL3 Polyclonal Antibody	A65070	Human	ELISA, WB, IHC
TACR1 Polyclonal Antibody	A54105	Human	ELISA, IHC, IF
TADA2B Polyclonal Antibody	A65754	Human	ELISA, IHC
TADA3 Polyclonal Antibody	A51933	Human	ELISA
TAF12 Polyclonal Antibody	A56964	Human	ELISA, IHC
TAF13 Polyclonal Antibody	A64690	Human	ELISA, IHC
TAGAP Polyclonal Antibody	A61226	Human	ELISA, IHC
TAGLN Polyclonal Antibody	A51476	Human	ELISA, WB, IHC
TAL1 Polyclonal Antibody	A68727	Human, Rat	ELISA, WB, IF
TALDO1 Polyclonal Antibody	A51941	Human	ELISA, WB, IHC
TANGO2 Polyclonal Antibody	A65422	Human	ELISA
TAPBPL Polyclonal Antibody	A67922	Human	ELISA, IHC
tar Polyclonal Antibody	A56867	Escherichia coli	ELISA, WB
TARS2 Polyclonal Antibody	A61230	Human	ELISA, IHC
TAS2R13 Polyclonal Antibody	A67998	Human	ELISA, IHC, IF
TAS2R14 Polyclonal Antibody	A67414	Human	ELISA, IHC
TAS2R41 Polyclonal Antibody	A63536	Human	ELISA, IHC, IF
TAS2R42 Polyclonal Antibody	A65742	Human	ELISA, IHC
TAX1BP1 Polyclonal Antibody	A68991	Human, Rat	ELISA, WB
TAX1BP3 Polyclonal Antibody	A51777	Human	ELISA, WB
TAZ Polyclonal Antibody	A63540	Human	ELISA, IHC, IF
TBC1D14 Polyclonal Antibody	A68010	Human	ELISA, WB, IHC
TBC1D19 Polyclonal Antibody	A61234	Human	ELISA, WB, IHC
TBC1D20 Polyclonal Antibody	A61238	Human, Mouse	ELISA, WB, IHC, IF
TBC1D22A Polyclonal Antibody	A61242	Human, Rat	ELISA, WB
TBC1D26 Polyclonal Antibody	A65806	Human, Mouse	ELISA, WB
TBC1D3 Polyclonal Antibody	A66130	Human	ELISA
TBC1D3C Polyclonal Antibody	A65378	Human	ELISA
TBC1D9B Polyclonal Antibody	A61246	Human	ELISA, IHC, IF
TBCA Polyclonal Antibody	A61250	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
TBCB Polyclonal Antibody	A61254	Human, Mouse	ELISA, WB, IHC, IP, IF
TBCC Polyclonal Antibody	A53617	Human, Mouse	ELISA, WB, IHC
TBCEL Polyclonal Antibody	A56555	Human, Rat	ELISA, WB
TBPL1 Polyclonal Antibody	A50516	Human	ELISA, IHC
TBX15 Polyclonal Antibody	A61258	Human	ELISA, WB
Tbx15 Polyclonal Antibody	A63544	Mouse	ELISA
TBX18 Polyclonal Antibody	A55599	Human	ELISA
TCEAL7 Polyclonal Antibody	A61262	Human	ELISA, WB, IHC
TCEAL8 Polyclonal Antibody	A66178	Human	ELISA, IHC
TCEAL9 Polyclonal Antibody	A67894	Human	ELISA
TCEANC2 Polyclonal Antibody	A51113	Human	ELISA, IHC
TCEB1 Polyclonal Antibody	A54228	Human	ELISA
TCHP Polyclonal Antibody	A61266	Human	ELISA, IHC
TCN1 Polyclonal Antibody	A61270	Human	ELISA, IHC
TCN2 Polyclonal Antibody	A63548	Human	ELISA, WB, IF
TCP10 Polyclonal Antibody	A61274	Human	ELISA
TCP10L Polyclonal Antibody	A66414	Human	ELISA, IHC
TCP11L2 Polyclonal Antibody	A61278	Human	ELISA, WB, IHC
TCTA Polyclonal Antibody	A63552	Human, Mouse	ELISA, WB, IHC
TCTE1 Polyclonal Antibody	A61282	Human	ELISA, WB, IHC, IF
TCTEX1D2 Polyclonal Antibody	A61286	Human	ELISA, IHC
tcyA Polyclonal Antibody	A64326	Bacillus subtilis	ELISA
TDGF1 Polyclonal Antibody	A54133	Human	ELISA
TDGF1 Polyclonal Antibody	A54137	Human	ELISA
TDRD9 Polyclonal Antibody	A68290	Human	ELISA, IHC
TEAD3 Polyclonal Antibody	A55467	Human, Mouse	ELISA, WB
TEAD4 Polyclonal Antibody	A56599	Human, Mouse	ELISA, WB, IHC
TEKT3 Polyclonal Antibody	A67450	Human	ELISA, IHC
TEKT4 Polyclonal Antibody	A61290	Human	ELISA
TEKT5 Polyclonal Antibody	A56988	Human, Mouse	ELISA, WB
TERT Polyclonal Antibody	A53505	Human	ELISA, IHC
TES-26 Polyclonal Antibody	A57029	Toxocara canis	ELISA
TESC Polyclonal Antibody	A57293	Human, Mouse	ELISA, WB
TESPA1 Polyclonal Antibody	A62818	Human	ELISA, IHC, IF
TET3 Polyclonal Antibody	A50520	Human	ELISA, IHC
tetR Polyclonal Antibody	A64538	Escherichia coli	ELISA
tetX Polyclonal Antibody	A50817	Clostridium tetani	ELISA
TEX12 Polyclonal Antibody	A53276	Human	ELISA
TEX13A Polyclonal Antibody	A67794	Human	ELISA, IHC
Tex35 Polyclonal Antibody	A65318	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
TEX43 Polyclonal Antibody	A65726	Human	ELISA, IHC
TF Polyclonal Antibody	A51885	Human	ELISA, WB
TFAM Polyclonal Antibody	A50524	Human	ELISA, WB, IHC
TFB1M Polyclonal Antibody	A69735	Human	ELISA, WB, IHC, IF
TFEC Polyclonal Antibody	A61294	Human	ELISA, WB
TFF3 Polyclonal Antibody	A55026	Human	ELISA
TFF3 Polyclonal Antibody	A55137	Human	ELISA
TFRC Polyclonal Antibody	A57862	Human	ELISA, WB, IHC, IF
TG Polyclonal Antibody	A64144	Human	ELISA
TGFA Polyclonal Antibody	A54053	Human	ELISA, WB, IHC, IF
TGFB1 Polyclonal Antibody	A53563	Human	ELISA
TGFB1 Polyclonal Antibody	A56911	Bovine	ELISA
TGFB1 Polyclonal Antibody	A56951	Chicken	ELISA
tgfb1 Polyclonal Antibody	A63556	Oncorhynchus mykiss, Human	ELISA, WB
TGFB2 Polyclonal Antibody	A52030	Human	ELISA, WB
TGFB2 Polyclonal Antibody	A56435	Human, Mouse	ELISA, WB, IHC
TGFB3 Polyclonal Antibody	A52344	Human	ELISA, IHC
TGFB3 Polyclonal Antibody	A61298	Human	ELISA, IHC
TGFBRAP1 Polyclonal Antibody	A66250	Human	ELISA, IHC
TGM2 Polyclonal Antibody	A55125	Human	ELISA, WB
TGM4 Polyclonal Antibody	A56543	Human	ELISA
TGS1 Polyclonal Antibody	A51145	Human, Mouse	ELISA, WB, IHC
THAP11 Polyclonal Antibody	A56943	Human	ELISA
THAP12 Polyclonal Antibody	A55551	Human	ELISA, WB, IHC
Thbs1 Polyclonal Antibody	A57733	Mouse	ELISA
THBS1 Polyclonal Antibody	A57801	Human	ELISA
Thbs2 Polyclonal Antibody	A57747	Mouse	ELISA
Thbs3 Polyclonal Antibody	A57726	Mouse	ELISA
Thbs3 Polyclonal Antibody	A57777	Mouse, Human	ELISA, WB
THEMIS Polyclonal Antibody	A69119	Human, Rat	ELISA, WB, IF
thiO Polyclonal Antibody	A56603	Bacillus subtilis	ELISA
THOC3 Polyclonal Antibody	A66302	Human	ELISA, IHC
THOC5 Polyclonal Antibody	A68823	Human	ELISA, IHC, IF
THOC7 Polyclonal Antibody	A65462	Human	ELISA
Thpo Polyclonal Antibody	A64108	Rat, Mouse	ELISA, WB
THRA Polyclonal Antibody	A50528	Human	ELISA, WB, IHC
THRB Polyclonal Antibody	A50532	Human, Mouse	ELISA, WB, IHC
THTPA Polyclonal Antibody	A61302	Human	ELISA, IHC
THY1 Polyclonal Antibody	A61306	Human	ELISA
TIFA Polyclonal Antibody	A61310	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
tig Polyclonal Antibody	A64570	Escherichia coli O127:H6	ELISA
TIGD4 Polyclonal Antibody	A66174	Human	ELISA, IHC
TIGD7 Polyclonal Antibody	A65606	Human	ELISA, IHC
TIGIT Polyclonal Antibody	A64998	Human	ELISA, IF
TIMELESS Polyclonal Antibody	A70339	Human	ELISA, IHC
TIMM10 Polyclonal Antibody	A50536	Human	ELISA, IHC
TIMM10B Polyclonal Antibody	A68162	Human, Mouse	ELISA, WB, IHC
TIMM8B Polyclonal Antibody	A61314	Human	ELISA, IHC
TIMM9 Polyclonal Antibody	A61318	Human	ELISA, IHC
Timp1 Polyclonal Antibody	A55379	Rat	ELISA
Timp1 Polyclonal Antibody	A63560	Mouse	ELISA
TIMP2 Polyclonal Antibody	A51769	Human	ELISA, WB
TIMP3 Polyclonal Antibody	A52050	Human	ELISA, WB
TIMP4 Polyclonal Antibody	A53017	Human	ELISA, IHC, IF
TIMP4 Polyclonal Antibody	A57089	Bovine	ELISA
TINAGL1 Polyclonal Antibody	A67558	Human	ELISA, IHC, IF
TIPRL Polyclonal Antibody	A61322	Human	ELISA
TK1 Polyclonal Antibody	A50540	Human	ELISA, WB, IHC
TK1 Polyclonal Antibody	A63564	Human herpesvirus 1	ELISA
TKFC Polyclonal Antibody	A68907	Human	ELISA, IF
TKTL2 Polyclonal Antibody	A67482	Human	ELISA, IHC
TLE4 Polyclonal Antibody	A68811	Human	ELISA, IF
TLN1 Polyclonal Antibody	A56463	Human	ELISA, IHC
TLN2 Polyclonal Antibody	A55639	Human	ELISA, WB, IHC
TLR2 Polyclonal Antibody	A54840	Human	ELISA
TLR6 Polyclonal Antibody	A61326	Human	ELISA, WB, IHC
Tlr7 Polyclonal Antibody	A53663	Mouse, Human	ELISA, WB
TM2D3 Polyclonal Antibody	A67222	Human	ELISA, IHC, IF
TM4SF1 Polyclonal Antibody	A61330	Human	ELISA, IHC, IF
TM4SF5 Polyclonal Antibody	A63568	Human	ELISA, IHC
TM7SF3 Polyclonal Antibody	A61334	Human	ELISA, IHC, IF
TMBIM1 Polyclonal Antibody	A61338	Human, Mouse	ELISA, WB, IF
TMCC1 Polyclonal Antibody	A63572	Human, Mouse	ELISA, WB, IHC, IF
TMCO5A Polyclonal Antibody	A61342	Human	ELISA, WB, IHC, IF
TMED2 Polyclonal Antibody	A61346	Human, Mouse	ELISA, WB, IHC
TMED3 Polyclonal Antibody	A61350	Human	ELISA, WB, IHC
TMED9 Polyclonal Antibody	A53113	Human, Mouse	ELISA, WB
TMEM100 Polyclonal Antibody	A61354	Human	ELISA, IF
TMEM106A Polyclonal Antibody	A61358	Human	ELISA, IF
TMEM109 Polyclonal Antibody	A64068	Human	ELISA, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
TMEM115 Polyclonal Antibody	A61362	Human	ELISA, IHC, IF
TMEM120B Polyclonal Antibody	A63576	Human	ELISA, IHC, IF
TMEM139 Polyclonal Antibody	A63580	Human	ELISA, IHC, IF
TMEM140 Polyclonal Antibody	A63584	Human	ELISA, IHC, IF
TMEM141 Polyclonal Antibody	A50544	Human	ELISA, WB, IHC
TMEM143 Polyclonal Antibody	A63588	Human	ELISA, IHC, IF
TMEM14A Polyclonal Antibody	A61366	Human	ELISA, WB, IF
TMEM14B Polyclonal Antibody	A57878	Human	ELISA, WB, IHC, IF
TMEM154 Polyclonal Antibody	A63592	Human	ELISA, WB, IHC
TMEM159 Polyclonal Antibody	A51157	Human	ELISA, IF
TMEM161A Polyclonal Antibody	A61370	Human	ELISA, WB, IHC, IF
TMEM161B Polyclonal Antibody	A63596	Human	ELISA
TMEM165 Polyclonal Antibody	A61374	Human	ELISA, WB, IHC, IF
TMEM174 Polyclonal Antibody	A61378	Human, Mouse	ELISA, WB, IHC, IF
TMEM176B Polyclonal Antibody	A61382	Human	ELISA, WB, IHC, IF
TMEM177 Polyclonal Antibody	A63600	Human	ELISA, WB, IHC, IF
TMEM184B Polyclonal Antibody	A61386	Human	ELISA, IHC, IF
TMEM184C Polyclonal Antibody	A63604	Human	ELISA, IHC, IF
TMEM186 Polyclonal Antibody	A61390	Human, Mouse	ELISA, WB, IHC, IF
TMEM19 Polyclonal Antibody	A63608	Human, Mouse	ELISA, WB, IHC, IF
TMEM192 Polyclonal Antibody	A61394	Human	ELISA, WB, IHC
TMEM196 Polyclonal Antibody	A63612	Human	ELISA
TMEM205 Polyclonal Antibody	A63616	Human	ELISA, IHC
TMEM207 Polyclonal Antibody	A63620	Human	ELISA, IHC, IF
TMEM209 Polyclonal Antibody	A63624	Human	ELISA, IHC, IF
TMEM213 Polyclonal Antibody	A63628	Human	ELISA, IHC, IF
TMEM217 Polyclonal Antibody	A61398	Human	ELISA, WB, IF
TMEM222 Polyclonal Antibody	A61402	Human	ELISA, WB, IHC, IF
TMEM243 Polyclonal Antibody	A67534	Human	ELISA, IHC, IF
TMEM25 Polyclonal Antibody	A63632	Human	ELISA
TMEM255A Polyclonal Antibody	A65130	Human, Mouse	ELISA, WB, IF
TMEM255B Polyclonal Antibody	A61406	Human	ELISA, IHC, IF
TMEM256 Polyclonal Antibody	A61410	Human	ELISA, IF
TMEM257 Polyclonal Antibody	A62354	Human	ELISA, IHC
TMEM27 Polyclonal Antibody	A51353	Human	ELISA, IHC, IF
TMEM31 Polyclonal Antibody	A63636	Human	ELISA, IHC, IF
TMEM33 Polyclonal Antibody	A61414	Human	ELISA, IHC, IF
TMEM38B Polyclonal Antibody	A61418	Human, Mouse	ELISA, WB, IHC, IF
TMEM40 Polyclonal Antibody	A63640	Human	ELISA, IHC, IF
TMEM41A Polyclonal Antibody	A61422	Human, Mouse	ELISA, WB, IHC, IF



Product Name	Cat. #	Species Reactivity	Applications
TMEM44 Polyclonal Antibody	A56875	Human	ELISA, IHC, IF
TMEM45A Polyclonal Antibody	A61426	Human	ELISA, IHC, IF
TMEM47 Polyclonal Antibody	A63644	Human	ELISA, IHC, IF
TMEM52B Polyclonal Antibody	A65034	Human	ELISA, IHC
TMEM53 Polyclonal Antibody	A63648	Human, Mouse	ELISA, WB, IHC, IF
TMEM55A Polyclonal Antibody	A61430	Human, Mouse	ELISA, WB, IHC, IF
TMEM56 Polyclonal Antibody	A63652	Human	ELISA, IHC, IF
TMEM59L Polyclonal Antibody	A61434	Human	ELISA, WB, IHC, IF
TMEM61 Polyclonal Antibody	A61438	Human	ELISA, IHC, IF
TMEM63A Polyclonal Antibody	A61442	Human	ELISA, IF
TMEM65 Polyclonal Antibody	A63656	Human, Mouse	ELISA, WB, IHC, IF
TMEM67 Polyclonal Antibody	A65090	Human	ELISA, IHC, IF
TMEM68 Polyclonal Antibody	A63660	Human, Mouse	ELISA, WB, IF
TMEM71 Polyclonal Antibody	A63664	Human	ELISA, WB, IHC
TMEM82 Polyclonal Antibody	A63668	Human	ELISA, IHC, IF
TMEM9 Polyclonal Antibody	A61446	Human	ELISA, IHC
TMEM91 Polyclonal Antibody	A63672	Human	ELISA, WB, IHC, IF
TMEM95 Polyclonal Antibody	A63676	Human	ELISA, IHC, IF
TMEM98 Polyclonal Antibody	A57373	Human	ELISA
tmk Polyclonal Antibody	A56579	Bacillus anthracis	ELISA
TMLHE Polyclonal Antibody	A54914	Human	ELISA, WB, IHC
TMPRSS11B Polyclonal Antibody	A65842	Human	ELISA, IHC
TMPRSS11D Polyclonal Antibody	A63680	Human	ELISA, IHC
Tmprss4 Polyclonal Antibody	A51121	Mouse	ELISA
TMPRSS6 Polyclonal Antibody	A61450	Human	ELISA, IHC
TMPRSS7 Polyclonal Antibody	A65458	Human	ELISA, IHC
TMSB4X Polyclonal Antibody	A69739	Human	ELISA, IHC
TMSB4Y Polyclonal Antibody	A63684	Human	ELISA
TMTC1 Polyclonal Antibody	A66014	Human, Mouse	ELISA, WB, IF
TMTC4 Polyclonal Antibody	A61454	Human	ELISA, IHC
TMX2 Polyclonal Antibody	A63688	Human	ELISA, IHC, IF
TMX3 Polyclonal Antibody	A55583	Human	ELISA, IHC
TNC Polyclonal Antibody	A52217	Human	ELISA, IHC
TNC Polyclonal Antibody	A52305	Human	ELISA, IHC
TNF Polyclonal Antibody	A53670	Sheep	ELISA, WB
TNF Polyclonal Antibody	A54089	Bovine	ELISA, WB
TNF Polyclonal Antibody	A54612	Human, Mouse	ELISA, WB, IHC
TNFAIP6 Polyclonal Antibody	A55182	Human	ELISA, IHC
TNFRSF10D Polyclonal Antibody	A54081	Human	ELISA, IHC
TNFRSF11A Polyclonal Antibody	A51841	Human	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
TNFRSF13C Polyclonal Antibody	A51181	Human	ELISA
TNFRSF19 Polyclonal Antibody	A69239	Human	ELISA, WB, IF
Tnfrsf4 Polyclonal Antibody	A61458	Mouse	ELISA
TNFSF10 Polyclonal Antibody	A51937	Human	ELISA, IHC
Tnfsf11 Polyclonal Antibody	A54073	Mouse	ELISA
Tnfsf11 Polyclonal Antibody	A63692	Mouse, Human	ELISA, IF
TNFSF12 Polyclonal Antibody	A61462	Human	ELISA, WB, IHC
TNFSF13B Polyclonal Antibody	A55647	Human	ELISA
TNFSF14 Polyclonal Antibody	A51761	Human	ELISA, WB, IF
TNFSF9 Polyclonal Antibody	A51974	Human	ELISA, IHC
TNIP1 Polyclonal Antibody	A68863	Human, Rat	ELISA, WB, IHC, IF
TNKS1BP1 Polyclonal Antibody	A69219	Human	ELISA, IHC, IF
TNNI3 Polyclonal Antibody	A51521	Human	ELISA
TNNI3 Polyclonal Antibody	A53028	Human	ELISA, IHC
Tnnt2 Polyclonal Antibody	A50548	Mouse	ELISA
TNNT2 Polyclonal Antibody	A52106	Human, Mouse, Rat	ELISA, WB, IHC
TNRC6A Polyclonal Antibody	A68282	Human	ELISA, IHC
TNRC6B Polyclonal Antibody	A69323	Human	ELISA, IF
TOX Polyclonal Antibody	A68731	Human	ELISA, WB
toxA Polyclonal Antibody	A64294	Peptoclostridium difficile	ELISA
toxB Polyclonal Antibody	A50760	Peptoclostridium difficile	ELISA
Toxin Zeta Polyclonal Antibody	A55034	Streptococcus agalactiae	ELISA
Tp53 Polyclonal Antibody	A52070	Rat, Mouse	ELISA, WB
Tp53 Polyclonal Antibody	A57385	Mouse	ELISA
TP53 Polyclonal Antibody	A68310	Human, Mouse	ELISA, WB, IHC, CHIP
TP53TG5 Polyclonal Antibody	A61466	Human	ELISA, IHC
Tpbg Polyclonal Antibody	A50552	Rat	ELISA
TPM2 Polyclonal Antibody	A53133	Human	ELISA
TPMT Polyclonal Antibody	A51617	Human	ELISA, WB, IHC
TPPP2 Polyclonal Antibody	A63696	Human, Mouse	ELISA, WB, IHC, IF
TPPP3 Polyclonal Antibody	A63700	Human	ELISA, IHC
TPRKB Polyclonal Antibody	A68174	Human	ELISA, WB, IHC
Tpsab1 Polyclonal Antibody	A53145	Mouse	ELISA
Tpsb2 Polyclonal Antibody	A53525	Mouse, Rat	ELISA, WB
Tpsb2 Polyclonal Antibody	A53574	Rat	ELISA
TPSB2 Polyclonal Antibody	A63704	Human, Mouse, Rat	ELISA, WB, IHC
TPST2 Polyclonal Antibody	A55229	Human	ELISA, IHC
Tpt1 Polyclonal Antibody	A52752	Mouse, Human	ELISA, IHC
TPTE Polyclonal Antibody	A50556	Human	ELISA, WB, IF
TPTE2 Polyclonal Antibody	A65626	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
TRAC Polyclonal Antibody	A55102	Human	ELISA
TRAF5 Polyclonal Antibody	A50560	Human	ELISA, IHC
TRAF6 Polyclonal Antibody	A54437	Human	ELISA, WB, IHC
TRAFD1 Polyclonal Antibody	A63708	Human	ELISA
TRAK1 Polyclonal Antibody	A68042	Human	ELISA, IHC
TRAM2 Polyclonal Antibody	A64846	Human	ELISA, IHC, IF
Transferrin Polyclonal Antibody	A57588	Rat	ELISA, WB
Transferrin Polyclonal Antibody	A57810	Sheep	ELISA
Transposase for Transposon Tn552 Polyclonal Antibody	A54332	Staphylococcus aureus	ELISA
TRAP1 Polyclonal Antibody	A51697	Human	ELISA, WB, IF
TRAPPC2L Polyclonal Antibody	A61470	Human, Mouse	ELISA, WB, IHC
TRAPPC3 Polyclonal Antibody	A61474	Human	ELISA, IHC
TRAPPC6A Polyclonal Antibody	A69743	Human	ELISA, IHC
TRAPPC6B Polyclonal Antibody	A65986	Human, Mouse	ELISA, WB, IHC
TRAT1 Polyclonal Antibody	A65706	Human	ELISA, IHC
TRBC1 Polyclonal Antibody	A56723	Human	ELISA, IHC, IF
Trem2 Polyclonal Antibody	A61478	Mouse, Human	ELISA, WB
TREML1 Polyclonal Antibody	A61482	Human	ELISA
TREML2 Polyclonal Antibody	A61486	Human	ELISA, IHC
TRHR Polyclonal Antibody	A68735	Human	ELISA, IF
TRIB1 Polyclonal Antibody	A69067	Human	ELISA, WB, IF
TRIM11 Polyclonal Antibody	A56122	Human	ELISA, IHC
TRIM15 Polyclonal Antibody	A61490	Human	ELISA, WB, IHC
TRIM16 Polyclonal Antibody	A63712	Human	ELISA
TRIM17 Polyclonal Antibody	A61494	Human	ELISA, WB, IHC
TRIM2 Polyclonal Antibody	A61498	Human	ELISA
Trim21 Polyclonal Antibody	A50985	Mouse	ELISA
TRIM28 Polyclonal Antibody	A51701	Human	ELISA, WB, IHC
TRIM35 Polyclonal Antibody	A68102	Human	ELISA
TRIM36 Polyclonal Antibody	A67398	Human	ELISA, IHC
TRIM37 Polyclonal Antibody	A63716	Human	ELISA, IHC, IF
TRIM38 Polyclonal Antibody	A61502	Human	ELISA, WB
TRIM43 Polyclonal Antibody	A66454	Human, Mouse	ELISA, WB
TRIM48 Polyclonal Antibody	A66162	Human	ELISA
TRIM51 Polyclonal Antibody	A61506	Human	ELISA, WB, IHC
TRIM55 Polyclonal Antibody	A67666	Human, Mouse	ELISA, WB, IHC
TRIM71 Polyclonal Antibody	A68899	Human	ELISA, IHC
TRIM74 Polyclonal Antibody	A61510	Human	ELISA, IHC
TRIM8 Polyclonal Antibody	A69215	Human	ELISA, IF
TRIM9 Polyclonal Antibody	A51345	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
TRIML1 Polyclonal Antibody	A61514	Human, Mouse	ELISA, WB
TRIP13 Polyclonal Antibody	A61518	Human	ELISA, WB, IHC
TRIT1 Polyclonal Antibody	A61522	Human	ELISA, WB, IHC
TRMT112 Polyclonal Antibody	A55451	Human	ELISA, IHC
TRMT6 Polyclonal Antibody	A68066	Human	ELISA, WB, IHC
TRNAU1AP Polyclonal Antibody	A61526	Human	ELISA, WB, IHC
TROVE2 Polyclonal Antibody	A52457	Human, Mouse, Rat	ELISA, WB, IHC, IF
TRPM2 Polyclonal Antibody	A62494	Human, Mouse	ELISA, WB, IHC, IF
TRPM8 Polyclonal Antibody	A51013	Human	ELISA, IHC, IF
TRPV2 Polyclonal Antibody	A69355	Human, Mouse	ELISA, WB, IF
TRPV6 Polyclonal Antibody	A70187	Human	ELISA, IHC
TRUB1 Polyclonal Antibody	A61530	Human	ELISA
trxA Polyclonal Antibody	A57734	Escherichia coli	ELISA, WB
trxC Polyclonal Antibody	A64258	Escherichia coli	ELISA
Trypsin Inhibitor DE-3 Polyclonal Antibody	A56919	Erythrina caffra	ELISA
Trypsin Inhibitor DE-3 Polyclonal Antibody	A64518	Erythrina caffra	ELISA
TSACC Polyclonal Antibody	A61534	Human	ELISA, IHC
TSC22D1 Polyclonal Antibody	A61538	Human	ELISA
TSC22D3 Polyclonal Antibody	A67158	Human	ELISA, WB, IHC
TSGA10 Polyclonal Antibody	A61542	Human	ELISA, WB, IHC
TSGA10IP Polyclonal Antibody	A64922	Human	ELISA
TSGA13 Polyclonal Antibody	A61546	Human	ELISA, WB, IHC
TSHB Polyclonal Antibody	A50684	Human	ELISA
TSHR Polyclonal Antibody	A53714	Human	ELISA, IF
TSLP Polyclonal Antibody	A63720	Human, Mouse	ELISA, WB, IHC
TSPAN1 Polyclonal Antibody	A61550	Human	ELISA, IHC, IF
TSPAN14 Polyclonal Antibody	A61554	Human	ELISA, IHC, IF
TSPAN15 Polyclonal Antibody	A61558	Human	ELISA, IF
TSPAN16 Polyclonal Antibody	A61562	Human	ELISA, IF
TSPAN31 Polyclonal Antibody	A50933	Human	ELISA, IHC, IF
TSPAN32 Polyclonal Antibody	A66726	Human	ELISA, IHC, IF
TSPAN33 Polyclonal Antibody	A65758	Human	ELISA, IHC, IF
TSPAN5 Polyclonal Antibody	A61566	Human	ELISA, WB, IF
TSPAN6 Polyclonal Antibody	A61570	Human, Mouse	ELISA, WB, IHC, IF
TSPAN7 Polyclonal Antibody	A54511	Human	ELISA, IHC, IF
TSPAN8 Polyclonal Antibody	A63724	Human, Mouse	ELISA, WB, IHC, IF
TSPAN9 Polyclonal Antibody	A68739	Human	ELISA, IHC, IF
TSSC1 Polyclonal Antibody	A61574	Human	ELISA, IHC
TST Polyclonal Antibody	A69887	Human, Rat	ELISA, WB, IHC
TSTA3 Polyclonal Antibody	A53877	Human	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
TTC1 Polyclonal Antibody	A52684	Human	ELISA, WB
TTC12 Polyclonal Antibody	A61578	Human, Mouse	ELISA, WB, IHC
TTC16 Polyclonal Antibody	A61582	Human	ELISA, IHC
TTC30A Polyclonal Antibody	A65994	Human, Mouse	ELISA, WB, IHC
TTC30B Polyclonal Antibody	A61586	Human	ELISA, WB, IHC
TTC32 Polyclonal Antibody	A65058	Human	ELISA, IHC
TTC33 Polyclonal Antibody	A61590	Human	ELISA, IHC
TTC38 Polyclonal Antibody	A61594	Human, Mouse	ELISA, WB, IHC
TTC5 Polyclonal Antibody	A66330	Human, Mouse	ELISA, WB
TTC9C Polyclonal Antibody	A66490	Human	ELISA, WB, IHC
TTF1 Polyclonal Antibody	A52101	Human	ELISA
TTL Polyclonal Antibody	A61598	Human	ELISA
TLL2 Polyclonal Antibody	A67226	Human	ELISA, IHC
TLL5 Polyclonal Antibody	A61602	Human	ELISA
TLL6 Polyclonal Antibody	A66198	Human	ELISA, IHC
TLL9 Polyclonal Antibody	A64962	Human	ELISA, IHC
TTN Polyclonal Antibody	A52472	Human	ELISA, IHC
Ttr Polyclonal Antibody	A53587	Mouse, Human	ELISA, IHC
TTR Polyclonal Antibody	A55707	Human	ELISA, IHC
Ttr Polyclonal Antibody	A56190	Mouse	ELISA
TTR Polyclonal Antibody	A57297	Macaca fascicularis	ELISA
TTYH1 Polyclonal Antibody	A61606	Human	ELISA, WB, IF
TUBA1A Polyclonal Antibody	A61610	Human	ELISA, IHC
TUBA3C Polyclonal Antibody	A64794	Human	ELISA
TUBA3E Polyclonal Antibody	A65582	Human	ELISA, WB, IHC
TUBAL3 Polyclonal Antibody	A63728	Human	ELISA, IHC
TUBB1 Polyclonal Antibody	A67370	Human, Mouse	ELISA, WB, IHC
TUBB2A Polyclonal Antibody	A56583	Human	ELISA, WB
TUBB3 Polyclonal Antibody	A51998	Human, Mouse	ELISA, WB, IHC
TUBB3 Polyclonal Antibody	A53965	Human, Mouse	ELISA, WB, IHC
TUBB6 Polyclonal Antibody	A51341	Human	ELISA, WB, IHC
TUBG2 Polyclonal Antibody	A61614	Human, Mouse	ELISA, WB, IHC
Tubulin Beta Chain A Polyclonal Antibody	A57753	Human	ELISA, IHC
Tubulin Beta Chain A Polyclonal Antibody	A57754	Human, Mouse, Zebrafish	ELISA, WB
Tubulin Beta-1 Chain Polyclonal Antibody	A52428	Gallus gallus	ELISA
tuf1 Polyclonal Antibody	A50564	Acinetobacter baumannii	ELISA
TUFM Polyclonal Antibody	A52852	Human	ELISA, WB
TULP3 Polyclonal Antibody	A63732	Human	ELISA, IHC
TWNK Polyclonal Antibody	A69095	Human, Mouse	ELISA, WB, IHC
TXN Polyclonal Antibody	A52015	Human	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
TXNDC11 Polyclonal Antibody	A61618	Human	ELISA, IHC, IF
TXNDC12 Polyclonal Antibody	A50568	Human	ELISA
TXNDC8 Polyclonal Antibody	A65194	Human	ELISA, IHC
TXNDC9 Polyclonal Antibody	A63736	Human	ELISA, IHC
Txnip Polyclonal Antibody	A66006	Mouse	ELISA
TXNIP Polyclonal Antibody	A69247	Human	ELISA, IF
TXNL1 Polyclonal Antibody	A54982	Human	ELISA, IHC
TXNRD1 Polyclonal Antibody	A50897	Human	ELISA, WB, IHC
TYMP Polyclonal Antibody	A63740	Human	ELISA, IHC
TYR Polyclonal Antibody	A56322	Human	ELISA, IHC
U2AF2 Polyclonal Antibody	A69747	Human	ELISA, WB
UBA5 Polyclonal Antibody	A61622	Human	ELISA, WB
UBA7 Polyclonal Antibody	A61626	Human	ELISA
UBAP1 Polyclonal Antibody	A67530	Human	ELISA, IHC
UBD Polyclonal Antibody	A50572	Human	ELISA, IHC, IF
UBE2C Polyclonal Antibody	A50576	Human	ELISA, WB, IHC
UBE2D1 Polyclonal Antibody	A50580	Human	ELISA, WB, IHC
UBE2D3 Polyclonal Antibody	A56803	Human	ELISA
UBE2DNL Polyclonal Antibody	A66066	Human	ELISA
UBE2G1 Polyclonal Antibody	A69751	Human	ELISA, IHC, IF
UBE2G2 Polyclonal Antibody	A61630	Human	ELISA, WB, IHC
UBE2I Polyclonal Antibody	A52612	Human	ELISA, WB, IHC
UBE2K Polyclonal Antibody	A53456	Human	ELISA, IHC
UBE2K Polyclonal Antibody	A57858	Human	ELISA
UBE2O Polyclonal Antibody	A61634	Human	ELISA, WB
UBE2Q2 Polyclonal Antibody	A54499	Human	ELISA, WB, IHC
UBE2T Polyclonal Antibody	A67594	Human	ELISA, WB, IHC
UBE2V1 Polyclonal Antibody	A63744	Human	ELISA, IHC
UBE2V2 Polyclonal Antibody	A55743	Human	ELISA, IHC
UBE2W Polyclonal Antibody	A55930	Human	ELISA
UBIAD1 Polyclonal Antibody	A69803	Human	ELISA, IF
UBL3 Polyclonal Antibody	A63748	Human	ELISA, IHC
UBL4A Polyclonal Antibody	A63752	Human	ELISA
UBL5 Polyclonal Antibody	A61638	Human	ELISA
UBL7 Polyclonal Antibody	A61642	Human	ELISA, IHC
UBLCP1 Polyclonal Antibody	A61646	Human	ELISA, WB, IHC
UBQLNL Polyclonal Antibody	A61650	Human	ELISA, IHC
UBR2 Polyclonal Antibody	A66030	Human, Mouse	ELISA, WB
UBTD1 Polyclonal Antibody	A61654	Human	ELISA
UBTD2 Polyclonal Antibody	A66790	Human	ELISA

Product Name	Cat. #	Species Reactivity	Applications
UBXN10 Polyclonal Antibody	A66470	Human	ELISA, IHC
UBXN2B Polyclonal Antibody	A64682	Human	ELISA, IHC, IF
UBXN6 Polyclonal Antibody	A61658	Human, Mouse	ELISA, WB, IHC
UBXN8 Polyclonal Antibody	A63756	Human	ELISA, IHC
UCHL1 Polyclonal Antibody	A52002	Human	ELISA
UCK1 Polyclonal Antibody	A55950	Human	ELISA, IHC
UCN2 Polyclonal Antibody	A66326	Human	ELISA, IHC
UCP1 Polyclonal Antibody	A54906	Human	ELISA, IHC
UCP2 Polyclonal Antibody	A63760	Human	ELISA
UCP3 Polyclonal Antibody	A63764	Human	ELISA, WB, IHC, IF
UFC1 Polyclonal Antibody	A61662	Human	ELISA
UFSP2 Polyclonal Antibody	A61666	Human, Mouse	ELISA, WB, IHC, IF
UGDH Polyclonal Antibody	A50584	Human	ELISA
UGP2 Polyclonal Antibody	A54856	Human, Mouse	ELISA, WB, IHC
UGT1A1 Polyclonal Antibody	A69755	Human, Mouse, Rat	ELISA, WB, IHC, IF
UGT1A6 Polyclonal Antibody	A69759	Human	ELISA, IHC
UGT2A3 Polyclonal Antibody	A65650	Human	ELISA, IHC
UGT2B4 Polyclonal Antibody	A61670	Human	ELISA, IHC
UHRF1BP1L Polyclonal Antibody	A61674	Human	ELISA
UL48 Polyclonal Antibody	A64442	Human herpesvirus 1	ELISA
UMPS Polyclonal Antibody	A68939	Human	ELISA, IF
UNC50 Polyclonal Antibody	A61678	Human	ELISA, WB, IF
UNC5B Polyclonal Antibody	A69763	Human	ELISA, IHC, IF
UNC93A Polyclonal Antibody	A65882	Human	ELISA, WB, IHC
UPP2 Polyclonal Antibody	A61682	Human, Mouse	ELISA, WB
UPRT Polyclonal Antibody	A61686	Human	ELISA, IHC
UQCR11 Polyclonal Antibody	A51501	Human	ELISA, IF
UQCERS1 Polyclonal Antibody	A51821	Human	ELISA, WB, IHC
UQCRQ Polyclonal Antibody	A51565	Human	ELISA, WB
ureA Polyclonal Antibody	A53606	Helicobacter pylori	ELISA, WB
ureB Polyclonal Antibody	A52062	Helicobacter pylori	ELISA, WB
UROS Polyclonal Antibody	A63768	Human	ELISA
USF1 Polyclonal Antibody	A54724	Human	ELISA, WB, IHC
USF2 Polyclonal Antibody	A68883	Human, Rat	ELISA, WB, IF
USH1C Polyclonal Antibody	A55850	Human	ELISA
USH1G Polyclonal Antibody	A65010	Human	ELISA
USHBP1 Polyclonal Antibody	A66634	Human	ELISA, IHC, IF
USMG5 Polyclonal Antibody	A61690	Human	ELISA, IF
USP12 Polyclonal Antibody	A50588	Human, Mouse	ELISA, WB
USP14 Polyclonal Antibody	A54547	Human	ELISA, WB, IHC, IF

Product Name	Cat. #	Species Reactivity	Applications
USP24 Polyclonal Antibody	A69303	Human	ELISA, IHC
USP3 Polyclonal Antibody	A68178	Human	ELISA, IHC
USP30 Polyclonal Antibody	A61694	Human, Mouse	ELISA, WB, IHC
USP33 Polyclonal Antibody	A66654	Human	ELISA, WB, IHC, IF
USP37 Polyclonal Antibody	A69007	Human	ELISA, IHC
USP47 Polyclonal Antibody	A61698	Human	ELISA, IHC
USP5 Polyclonal Antibody	A63772	Human, Mouse	ELISA, WB, IHC
uspF Polyclonal Antibody	A64378	Escherichia coli	ELISA
UTP15 Polyclonal Antibody	A61702	Human	ELISA, WB
UTP23 Polyclonal Antibody	A61706	Human, Mouse	ELISA, WB, IHC
UTP6 Polyclonal Antibody	A51289	Human	ELISA
UTY Polyclonal Antibody	A68743	Human	ELISA, IF
uvyY Polyclonal Antibody	A64542	Enterobacteria phage T4	ELISA
UXS1 Polyclonal Antibody	A61710	Human, Mouse	ELISA, WB
vacA Polyclonal Antibody	A52468	Helicobacter pylori	ELISA
vacA Polyclonal Antibody	A55020	Helicobacter pylori	ELISA
vacA Polyclonal Antibody	A64398	Helicobacter pylori	ELISA
VAMP1 Polyclonal Antibody	A63776	Human	ELISA, IHC
VAMP5 Polyclonal Antibody	A63780	Human	ELISA, IHC
VAMP7 Polyclonal Antibody	A52580	Chicken, Mouse	ELISA, IHC
VAMP7 Polyclonal Antibody	A54943	Human, Mouse	ELISA, WB, IHC, IF
VASH1 Polyclonal Antibody	A57185	Human	ELISA
VASH2 Polyclonal Antibody	A51017	Human	ELISA, IHC
VAT1L Polyclonal Antibody	A61714	Human	ELISA, WB, IHC
VAX1 Polyclonal Antibody	A69767	Human	ELISA, WB
VCAM1 Polyclonal Antibody	A55922	Human	ELISA
VCAM1 Polyclonal Antibody	A57890	Human	ELISA, IHC
VCL Polyclonal Antibody	A57633	Human	ELISA, WB
Vegfa Polyclonal Antibody	A52190	Rat, Mouse	ELISA, WB, IHC
VEGFA Polyclonal Antibody	A52744	Pig	ELISA
VEGFA Polyclonal Antibody	A61718	Human	ELISA, WB
VEGFC Polyclonal Antibody	A51773	Human	ELISA, WB, IHC
Vegfc Polyclonal Antibody	A54057	Mouse	ELISA
Vegfc Polyclonal Antibody	A57209	Mouse	ELISA
VGLL2 Polyclonal Antibody	A66498	Human	ELISA, IHC
VHLL Polyclonal Antibody	A65646	Human	ELISA, IHC
viaA Polyclonal Antibody	A53229	Salmonella typhi	ELISA
VIM Polyclonal Antibody	A63784	Human	ELISA, WB, IHC, IF
VIPAS39 Polyclonal Antibody	A61722	Human, Mouse	ELISA, WB
virD2 Polyclonal Antibody	A53969	Agrobacterium tumefaciens	ELISA



Product Name	Cat. #	Species Reactivity	Applications
virE2 Polyclonal Antibody	A64462	Agrobacterium fabrum	ELISA
VKORC1L1 Polyclonal Antibody	A66186	Human	ELISA, IHC
VMO1 Polyclonal Antibody	A65950	Human	ELISA, IHC, IF
VN1R2 Polyclonal Antibody	A66866	Human	ELISA, WB, IF
VN1R5 Polyclonal Antibody	A65978	Human	ELISA, IHC, IF
VOPP1 Polyclonal Antibody	A51733	Human	ELISA, WB
VOPP1 Polyclonal Antibody	A51737	Human	ELISA, WB, IF
VP35 Polyclonal Antibody	A67074	Reston ebolavirus	ELISA
VP40 Polyclonal Antibody	A66074	Reston ebolavirus	ELISA
VPREB3 Polyclonal Antibody	A61726	Human	ELISA
VPS25 Polyclonal Antibody	A67906	Human, Mouse	ELISA, WB, IHC
VPS26B Polyclonal Antibody	A61730	Human, Mouse	ELISA, WB, IHC, IF
VPS45 Polyclonal Antibody	A67730	Human, Mouse	ELISA, WB, IHC
VPS50 Polyclonal Antibody	A67106	Human	ELISA, IHC
VPS52 Polyclonal Antibody	A70147	Human	ELISA, IHC
vpx Polyclonal Antibody	A57053	Human immunodeficiency virus type 2 subtype A	ELISA
vpx Polyclonal Antibody	A64278	Simian immunodeficiency virus	ELISA
vpx Polyclonal Antibody	A64290	Simian immunodeficiency virus	ELISA
VSIG2 Polyclonal Antibody	A51097	Human	ELISA, IHC
VTA1 Polyclonal Antibody	A61734	Human	ELISA, IHC
VTCN1 Polyclonal Antibody	A52318	Human	ELISA
VTG2 Polyclonal Antibody	A53468	Gallus gallus	ELISA
VTI1B Polyclonal Antibody	A56254	Human	ELISA
VWA3B Polyclonal Antibody	A61738	Human	ELISA, IHC
VWA5A Polyclonal Antibody	A61742	Human	ELISA, WB
VWA5B2 Polyclonal Antibody	A61746	Human	ELISA, IHC
VWA9 Polyclonal Antibody	A66834	Human, Mouse	ELISA, WB, IHC
VWF Polyclonal Antibody	A55683	Human	ELISA, IHC
WASF2 Polyclonal Antibody	A69335	Human	ELISA, IHC, IF
WBP2 Polyclonal Antibody	A61750	Human	ELISA, IHC
WBSCR17 Polyclonal Antibody	A65426	Human	ELISA, IHC
WBSCR28 Polyclonal Antibody	A61754	Human	ELISA, IHC, IF
WDFY3 Polyclonal Antibody	A69995	Human	ELISA, IHC, IF
Wdpcp Polyclonal Antibody	A53333	Mouse	ELISA, WB
WDR18 Polyclonal Antibody	A61758	Human	ELISA
WDR27 Polyclonal Antibody	A63792	Human	ELISA, IHC
WDR34 Polyclonal Antibody	A66702	Human	ELISA, WB, IHC
Wdr37 Polyclonal Antibody	A69039	Human, Mouse	ELISA, WB

Product Name	Cat. #	Species Reactivity	Applications
WDR37 Polyclonal Antibody	A61762	Human	ELISA, WB, IHC
WDR49 Polyclonal Antibody	A61766	Human	ELISA, WB, IHC
WDR5 Polyclonal Antibody	A56831	Human	ELISA
WDR53 Polyclonal Antibody	A65858	Human	ELISA, IHC
WDR54 Polyclonal Antibody	A55315	Human, Mouse	ELISA, WB, IHC
WDR61 Polyclonal Antibody	A61770	Human, Mouse	ELISA, WB, IHC
WDR62 Polyclonal Antibody	A63796	Human	ELISA, IHC
WDR66 Polyclonal Antibody	A66874	Human	ELISA, IHC
WDR74 Polyclonal Antibody	A65486	Human	ELISA, WB, IHC
WDR77 Polyclonal Antibody	A56947	Human	ELISA, IHC
WDR88 Polyclonal Antibody	A61774	Human	ELISA, WB, IHC
WDR89 Polyclonal Antibody	A61778	Human	ELISA, IHC
WDR91 Polyclonal Antibody	A63800	Human	ELISA, IHC
WDTC1 Polyclonal Antibody	A61782	Human, Mouse	ELISA, WB
WDYHV1 Polyclonal Antibody	A61786	Human, Mouse	ELISA, WB
WFDC11 Polyclonal Antibody	A66222	Human	ELISA, IHC
WFDC2 Polyclonal Antibody	A51925	Human, Mouse	ELISA, WB, IHC, IF
WFDC9 Polyclonal Antibody	A66402	Human	ELISA
Wheat Gliadin Polyclonal Antibody	A57785	Plant	ELISA, WB
WHSC1L1 Polyclonal Antibody	A61790	Human	ELISA, IHC
WLS Polyclonal Antibody	A68919	Human	ELISA, WB, IHC
Wnt3 Polyclonal Antibody	A61794	Mouse, Human	ELISA, WB
WNT3A Polyclonal Antibody	A52869	Human, Mouse	ELISA, IHC
WNT4 Polyclonal Antibody	A61798	Human	ELISA, WB, IHC
wnt8 Polyclonal Antibody	A50800	Xenopus laevis	ELISA
WNT8A Polyclonal Antibody	A70207	Human	ELISA, IHC
WNT9B Polyclonal Antibody	A69775	Human	ELISA, IHC
WRKY40 Polyclonal Antibody	A54352	Arabidopsis thaliana	ELISA
XAGE1A Polyclonal Antibody	A61802	Human	ELISA, IHC
XAGE2 Polyclonal Antibody	A66594	Human	ELISA
XG Polyclonal Antibody	A63804	Human, Mouse	ELISA, WB
XIAP Polyclonal Antibody	A50592	Human	ELISA, IHC
XIRP2 Polyclonal Antibody	A69779	Human	ELISA, IHC
XPC Polyclonal Antibody	A53796	Human	ELISA
XPNPEP1 Polyclonal Antibody	A55603	Human, Mouse	ELISA, WB, IHC
XPR1 Polyclonal Antibody	A67426	Human	ELISA, IHC, IF
XRCC3 Polyclonal Antibody	A57289	Human	ELISA
XRCC5 Polyclonal Antibody	A61806	Human	ELISA, WB, IHC
XRCC6 Polyclonal Antibody	A56647	Human, Mouse	ELISA, WB, IHC, IF
XRRA1 Polyclonal Antibody	A65478	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
YAF2 Polyclonal Antibody	A66070	Human	ELISA, IHC
YARS Polyclonal Antibody	A52632	Human	ELISA, WB, IHC
ydhF Polyclonal Antibody	A53064	Escherichia coli	ELISA
yedQ Polyclonal Antibody	A51801	Escherichia coli	ELISA, WB
YIF1B Polyclonal Antibody	A50965	Human, Rat	ELISA, WB, IHC, IF
YIPF4 Polyclonal Antibody	A61810	Human	ELISA, WB, IF
YIPF5 Polyclonal Antibody	A61814	Human, Mouse	ELISA, WB, IF
YIPF6 Polyclonal Antibody	A61818	Human	ELISA, IHC, IF
YKT6 Polyclonal Antibody	A61822	Human	ELISA, WB, IHC
yoaJ Polyclonal Antibody	A57823	Bacillus subtilis	ELISA
yoaJ Polyclonal Antibody	A57838	Bacillus subtilis	ELISA
YPEL1 Polyclonal Antibody	A50596	Human	ELISA, IHC
YPEL3 Polyclonal Antibody	A57321	Human	ELISA, IHC
YPEL5 Polyclonal Antibody	A68747	Human	ELISA, IHC
YTHDF2 Polyclonal Antibody	A68190	Human	ELISA, WB, IHC, IF
YTHDF3 Polyclonal Antibody	A65954	Human	ELISA
YWHAB Polyclonal Antibody	A56655	Human, Mouse, Rat, Zebrafish	ELISA, WB, IHC, IF
YWHAE Polyclonal Antibody	A50600	Human, Mouse	ELISA, WB, IHC
YWHAH Polyclonal Antibody	A53158	Human, Mouse	ELISA, WB, IHC
ZAR1 Polyclonal Antibody	A68903	Arabidopsis thaliana	ELISA
ZBED1 Polyclonal Antibody	A53917	Human	ELISA, WB, IHC
ZBED6CL Polyclonal Antibody	A61826	Human	ELISA, IHC
ZBTB18 Polyclonal Antibody	A61830	Human	ELISA, WB, IHC
ZBTB32 Polyclonal Antibody	A61834	Human, Mouse	ELISA, WB, IHC
ZBTB43 Polyclonal Antibody	A61838	Human, Mouse	ELISA, WB, IHC
ZBTB6 Polyclonal Antibody	A61842	Human	ELISA, WB, IHC
ZBTB8OS Polyclonal Antibody	A66150	Human	ELISA, WB, IHC, IF
ZC3H10 Polyclonal Antibody	A61846	Human	ELISA
ZC3H12A Polyclonal Antibody	A61850	Human	ELISA, IHC
ZC3H7A Polyclonal Antibody	A61854	Human	ELISA, IHC
ZCCHC13 Polyclonal Antibody	A66562	Human	ELISA, IHC
ZCCHC17 Polyclonal Antibody	A61858	Human	ELISA
ZCCHC4 Polyclonal Antibody	A61862	Human	ELISA, WB, IHC
ZDHHC12 Polyclonal Antibody	A66298	Human	ELISA, WB, IHC, IF
ZDHHC15 Polyclonal Antibody	A66718	Human	ELISA, IHC, IF
ZDHHC2 Polyclonal Antibody	A61870	Human	ELISA, IHC, IF
ZDHHC20 Polyclonal Antibody	A65238	Human	ELISA, IF
ZDHHC23 Polyclonal Antibody	A61866	Human, Mouse	ELISA, WB, IHC, IF
ZDHHC3 Polyclonal Antibody	A61874	Human, Mouse	ELISA, WB, IHC, IF
ZDHHC4 Polyclonal Antibody	A61878	Human, Mouse	ELISA, WB, IHC

Product Name	Cat. #	Species Reactivity	Applications
ZDHC5 Polyclonal Antibody	A61882	Human	ELISA, IHC, IF
ZDHC6 Polyclonal Antibody	A61886	Human, Mouse	ELISA, WB, IHC, IF
ZFAND2A Polyclonal Antibody	A61890	Human	ELISA, IHC
ZFP36 Polyclonal Antibody	A68751	Human	ELISA, IF
ZFP64 Polyclonal Antibody	A61894	Human, Mouse	ELISA, WB
ZFP69B Polyclonal Antibody	A63996	Human	ELISA, IHC
ZFP82 Polyclonal Antibody	A66738	Human	ELISA
ZFX Polyclonal Antibody	A69783	Human	ELISA, IHC
ZG16 Polyclonal Antibody	A61898	Human	ELISA, IHC
ZGRF1 Polyclonal Antibody	A66002	Human	ELISA, IF
ZHX1-C8orf76 Polyclonal Antibody	A61902	Human, Mouse	ELISA, WB, IHC
ZIC5 Polyclonal Antibody	A69115	Human	ELISA, IHC
ZKSCAN2 Polyclonal Antibody	A65306	Human	ELISA, IHC
ZKSCAN4 Polyclonal Antibody	A61906	Human	ELISA
ZKSCAN5 Polyclonal Antibody	A70343	Human	ELISA, IHC
ZKSCAN7 Polyclonal Antibody	A61910	Human	ELISA
ZKSCAN8 Polyclonal Antibody	A61914	Human	ELISA
ZMAT2 Polyclonal Antibody	A52269	Human	ELISA
ZMAT4 Polyclonal Antibody	A61918	Human	ELISA, IHC
ZMAT5 Polyclonal Antibody	A51265	Human	ELISA, WB, IHC
ZMYM3 Polyclonal Antibody	A64678	Human	ELISA, IF
ZMYND19 Polyclonal Antibody	A61922	Human	ELISA, IHC
ZNF114 Polyclonal Antibody	A55611	Human	ELISA, IHC
ZNF131 Polyclonal Antibody	A63808	Human	ELISA, IHC
ZNF133 Polyclonal Antibody	A61926	Human	ELISA, WB, IHC, IF
ZNF155 Polyclonal Antibody	A57357	Human	ELISA, IHC
ZNF165 Polyclonal Antibody	A50604	Human	ELISA, IHC
ZNF175 Polyclonal Antibody	A55311	Human	ELISA, IHC
ZNF177 Polyclonal Antibody	A63812	Human	ELISA, IHC
ZNF184 Polyclonal Antibody	A61930	Human	ELISA, IHC
ZNF19 Polyclonal Antibody	A61934	Human, Mouse	ELISA, WB, IHC
ZNF212 Polyclonal Antibody	A50608	Human, Mouse	ELISA, WB, IHC
ZNF215 Polyclonal Antibody	A63816	Human	ELISA, WB, IHC
ZNF221 Polyclonal Antibody	A63820	Human	ELISA, IHC
ZNF223 Polyclonal Antibody	A63824	Human	ELISA
ZNF226 Polyclonal Antibody	A63828	Human	ELISA, IHC
ZNF227 Polyclonal Antibody	A63832	Human	ELISA
ZNF230 Polyclonal Antibody	A63836	Human	ELISA, IHC
ZNF232 Polyclonal Antibody	A61938	Human, Mouse	ELISA, WB, IHC
ZNF252P-AS1 Polyclonal Antibody	A64622	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
ZNF26 Polyclonal Antibody	A50612	Human	ELISA, IHC
ZNF260 Polyclonal Antibody	A63840	Human	ELISA, IHC
ZNF266 Polyclonal Antibody	A61942	Human	ELISA, IHC
ZNF274 Polyclonal Antibody	A66918	Human	ELISA, IHC
ZNF281 Polyclonal Antibody	A63844	Human	ELISA, IHC
ZNF285 Polyclonal Antibody	A56767	Human	ELISA
ZNF287 Polyclonal Antibody	A63848	Human	ELISA, IHC
ZNF300 Polyclonal Antibody	A63852	Human	ELISA
ZNF32 Polyclonal Antibody	A63856	Human	ELISA
ZNF329 Polyclonal Antibody	A63860	Human	ELISA
ZNF333 Polyclonal Antibody	A63864	Human	ELISA
ZNF366 Polyclonal Antibody	A63868	Human	ELISA, IHC, IF
ZNF384 Polyclonal Antibody	A53949	Human	ELISA, WB, IHC
ZNF385C Polyclonal Antibody	A63872	Human	ELISA
ZNF396 Polyclonal Antibody	A63876	Human	ELISA, IHC
ZNF397 Polyclonal Antibody	A61946	Human	ELISA, IHC
ZNF398 Polyclonal Antibody	A63880	Human	ELISA, WB
ZNF404 Polyclonal Antibody	A63884	Human	ELISA, IHC
ZNF410 Polyclonal Antibody	A63888	Human, Mouse	ELISA, WB
ZNF432 Polyclonal Antibody	A61950	Human	ELISA, IHC
ZNF433 Polyclonal Antibody	A63892	Human	ELISA
ZNF440 Polyclonal Antibody	A63896	Human	ELISA
ZNF454 Polyclonal Antibody	A63900	Human	ELISA, IHC
ZNF462 Polyclonal Antibody	A63904	Human	ELISA, IHC
ZNF484 Polyclonal Antibody	A63908	Human	ELISA, IHC
ZNF486 Polyclonal Antibody	A63912	Human	ELISA
ZNF488 Polyclonal Antibody	A63916	Human	ELISA, WB, IHC
ZNF490 Polyclonal Antibody	A61954	Human	ELISA, IHC
ZNF502 Polyclonal Antibody	A61958	Human	ELISA, IHC
ZNF513 Polyclonal Antibody	A63920	Human	ELISA, IHC
ZNF529 Polyclonal Antibody	A63924	Human	ELISA, WB, IHC, IF
ZNF530 Polyclonal Antibody	A63928	Human	ELISA
ZNF546 Polyclonal Antibody	A63932	Human	ELISA, WB
ZNF554 Polyclonal Antibody	A63936	Human	ELISA, WB
ZNF555 Polyclonal Antibody	A63940	Human, Mouse	ELISA, WB, IHC
ZNF559 Polyclonal Antibody	A61962	Human	ELISA, IHC
ZNF560 Polyclonal Antibody	A63944	Human	ELISA, IHC
ZNF562 Polyclonal Antibody	A63948	Human	ELISA, WB, IHC
ZNF563 Polyclonal Antibody	A63952	Human	ELISA, WB, IHC
ZNF564 Polyclonal Antibody	A63956	Human	ELISA, IHC

Product Name	Cat. #	Species Reactivity	Applications
ZNF582 Polyclonal Antibody	A63960	Human	ELISA, IHC
ZNF583 Polyclonal Antibody	A63964	Human	ELISA, IHC
ZNF586 Polyclonal Antibody	A63968	Human	ELISA, IHC
ZNF592 Polyclonal Antibody	A52006	Human	ELISA
ZNF595 Polyclonal Antibody	A61966	Human	ELISA
ZNF596 Polyclonal Antibody	A63972	Human	ELISA, IHC, IF
ZNF606 Polyclonal Antibody	A63976	Human	ELISA
ZNF610 Polyclonal Antibody	A61970	Human	ELISA, WB
ZNF614 Polyclonal Antibody	A61974	Human	ELISA, WB, IHC
ZNF621 Polyclonal Antibody	A63980	Human	ELISA
ZNF622 Polyclonal Antibody	A63984	Human	ELISA, IHC
ZNF624 Polyclonal Antibody	A63988	Human	ELISA, IHC, IF
ZNF625 Polyclonal Antibody	A61978	Human, Mouse	ELISA, WB
ZNF626 Polyclonal Antibody	A61982	Human	ELISA, WB, IHC
ZNF627 Polyclonal Antibody	A63992	Human	ELISA, IHC
ZNF662 Polyclonal Antibody	A64000	Human	ELISA, IHC
ZNF670 Polyclonal Antibody	A61986	Human	ELISA, WB, IHC
ZNF679 Polyclonal Antibody	A61990	Human	ELISA, IHC
ZNF680 Polyclonal Antibody	A64004	Human	ELISA
ZNF688 Polyclonal Antibody	A61994	Human	ELISA, IHC
ZNF689 Polyclonal Antibody	A64008	Human	ELISA, WB
ZNF699 Polyclonal Antibody	A61998	Human	ELISA, WB, IHC
ZNF703 Polyclonal Antibody	A69787	Human, Rat	ELISA, WB, IHC, IF
ZNF707 Polyclonal Antibody	A62002	Human, Mouse	ELISA, WB, IHC
ZNF71 Polyclonal Antibody	A67854	Human	ELISA, IHC
ZNF738 Polyclonal Antibody	A62006	Human	ELISA, IHC
ZNF75A Polyclonal Antibody	A64012	Human, Mouse	ELISA, WB
ZNF765 Polyclonal Antibody	A64016	Human	ELISA
ZNF767P Polyclonal Antibody	A65410	Human	ELISA, IHC
ZNF77 Polyclonal Antibody	A64020	Human	ELISA, IHC
ZNF774 Polyclonal Antibody	A64024	Human	ELISA, IHC
ZNF791 Polyclonal Antibody	A64028	Human	ELISA, IHC
ZNF800 Polyclonal Antibody	A64032	Human	ELISA, IHC
ZNF821 Polyclonal Antibody	A64036	Human	ELISA, WB, IHC
ZNF829 Polyclonal Antibody	A64040	Human	ELISA, WB
ZNF844 Polyclonal Antibody	A64044	Human	ELISA, IHC
ZNF91 Polyclonal Antibody	A53496	Human	ELISA, IHC
ZNF98 Polyclonal Antibody	A64048	Human	ELISA, IF
ZNRD1 Polyclonal Antibody	A56739	Human	ELISA, WB
ZNRF3 Polyclonal Antibody	A68038	Human	ELISA, IF

Product Name	Cat. #	Species Reactivity	Applications
ZBPB Polyclonal Antibody	A62010	Human	ELISA
ZRANB2 Polyclonal Antibody	A54601	Human	ELISA, IHC
ZSCAN16 Polyclonal Antibody	A62014	Human	ELISA
ZSCAN20 Polyclonal Antibody	A56002	Human	ELISA, IHC
ZSCAN21 Polyclonal Antibody	A52058	Human	ELISA
ZSCAN29 Polyclonal Antibody	A66034	Human	ELISA, IHC
ZSCAN32 Polyclonal Antibody	A62018	Human	ELISA, WB
ZSCAN9 Polyclonal Antibody	A62022	Human, Mouse	ELISA, WB, IHC
ZXDC Polyclonal Antibody	A64870	Human	ELISA