

HDAC1 Polyclonal Antibody

(Catalog # A73224)

Background

Histone acetylation and deacetylation, catalyzed by multisubunit complexes, play a key role in the regulation of eukaryotic gene expression. The protein encoded by this gene belongs to the histone deacetylase/acuc/apha family and is a component of the histone deacetylase complex. It also interacts with retinoblastoma tumor-suppressor protein and this complex is a key element in the control of cell proliferation and differentiation. Together with metastasis-associated protein-2, it deacetylates p53 and modulates its effect on cell growth and apoptosis.

Description

HDAC1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q13547

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 393-482 of human HDAC1 (NP_004955.2).

Storage

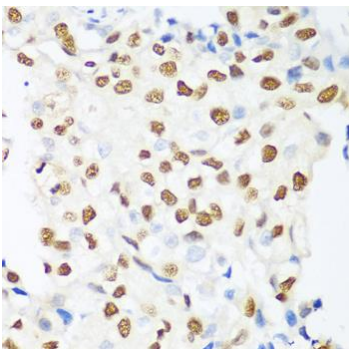
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

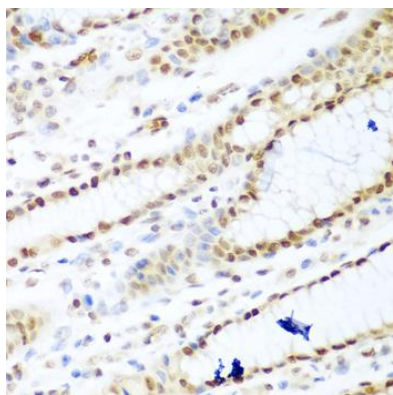
GON-10; HD1; RPD3; RPD3L1; HDAC1

Application

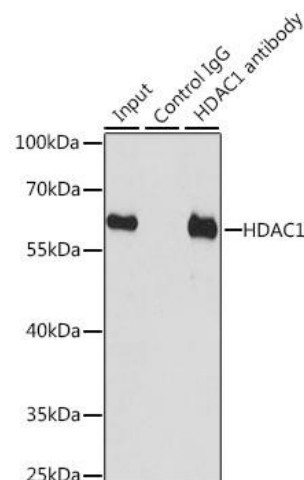
WB, IHC, IF, IP; Recommended dilution: WB 1:500 - 1:2000, IHC 1:100 - 1:200, IF 1:50 - 1:100, IP 1:50 - 1:100



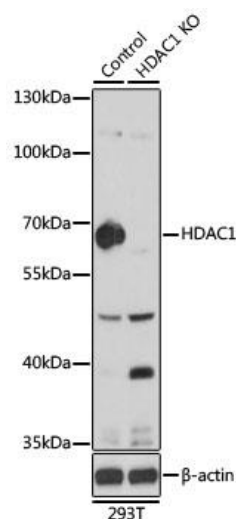
Immunohistochemistry of paraffin-embedded human prostate cancer using HDAC1 Polyclonal Antibody at dilution of 1:100 (40x lens).



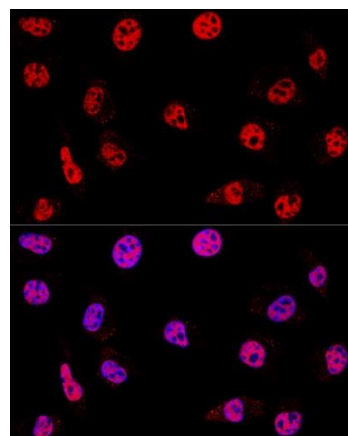
Immunohistochemistry of paraffin-embedded human stomach using HDAC1 Polyclonal Antibody at dilution of 1:100 (40x lens).



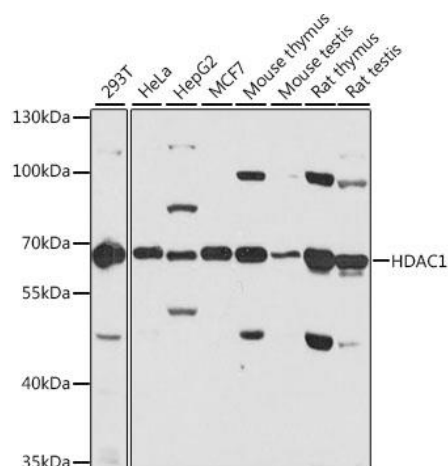
Immunoprecipitation analysis of 200ug extracts of HT-29 cells, using 3 ug HDAC1 Polyclonal Antibody. Western blot was performed from the immunoprecipitate using HDAC1 Polyclonal Antibody at a dilution of 1:1000.



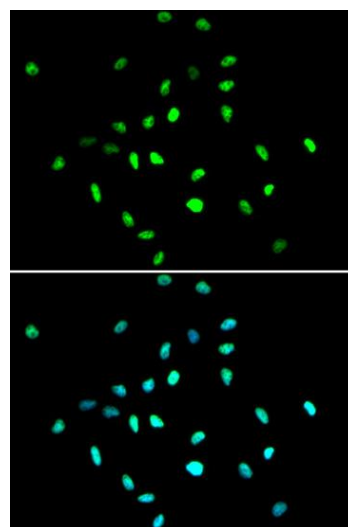
Western blot analysis of extracts from normal (control) and HDAC1 knockout (KO) 293T cells, using HDAC1 Polyclonal Antibody at 1:1000 dilution.
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 60s.



Confocal immunofluorescence analysis of L929 cells using HDAC1 Polyclonal Antibody at dilution of 1:200. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using HDAC1 Polyclonal Antibody at 1:1000 dilution.
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 60s.



Immunofluorescence analysis of A549 cells using HDAC1 Polyclonal Antibody. Blue: DAPI for nuclear staining.