
TRAF2 Polyclonal Antibody

(Catalog # A73233)

Background

The protein encoded by this gene is a member of the TNF receptor-associated factor (TRAF) protein family. TRAF proteins associate with and mediate the signal transduction from members of the TNF receptor superfamily. This protein directly interacts with TNF receptors and forms a heterodimeric complex with TRAF1. This protein is required for TNF-alpha-mediated activation of MAPK8/JNK and NF-kappaB. The protein complex formed by this protein and TRAF1 interacts with the inhibitor-of-apoptosis proteins (IAPs), and functions as a mediator of the anti-apoptotic signals from TNF receptors. The interaction of this protein with TRADD, a TNF receptor-associated apoptotic signal transducer, ensures the recruitment of IAPs for the direct inhibition of caspase activation. BIRC2/c-IAP1, an apoptosis inhibitor possessing ubiquitin ligase activity, can ubiquitinate and induce the degradation of this protein and thus potentiate TNF-induced apoptosis. Multiple alternatively spliced transcript variants have been found for this gene, but the biological validity of only one transcript has been determined.

Description

TRAF2 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

Q12933

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-300 of human TRAF2 (NP_066961.2).

Storage

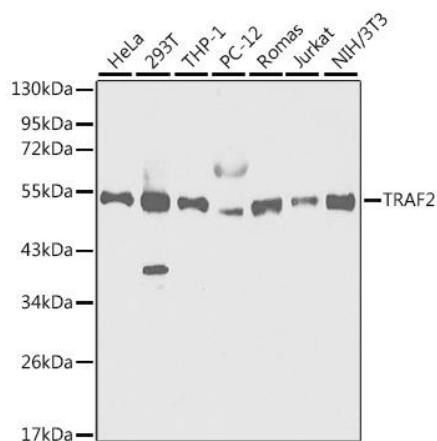
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

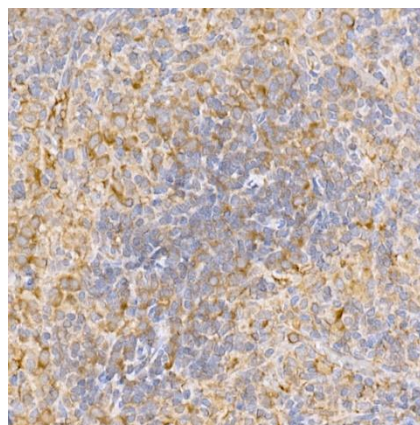
MGC:45012; TRAP; TRAP3; TRAF2

Application

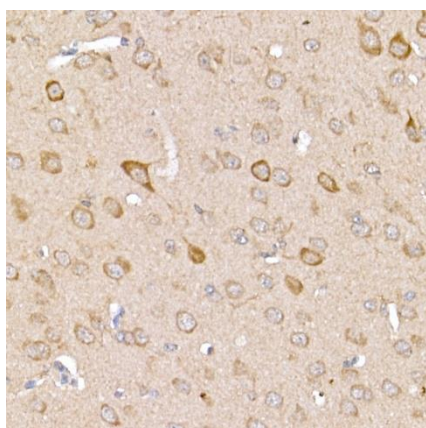
WB, IHC, IP; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:200, IP 1:50 - 1:100



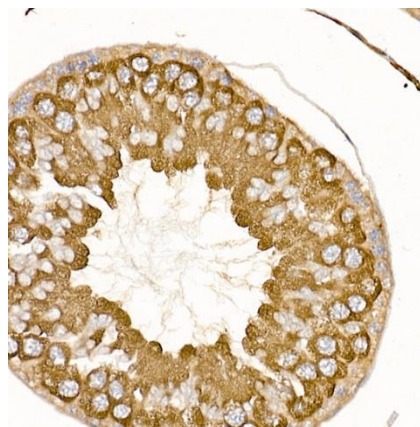
Western blot analysis of extracts of various cell lines, using TRAF2 Polyclonal Antibody at 1:1000 dilution.
 Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
 Lysates/proteins: 25ug per lane.
 Blocking buffer: 3% nonfat dry milk in TBST.



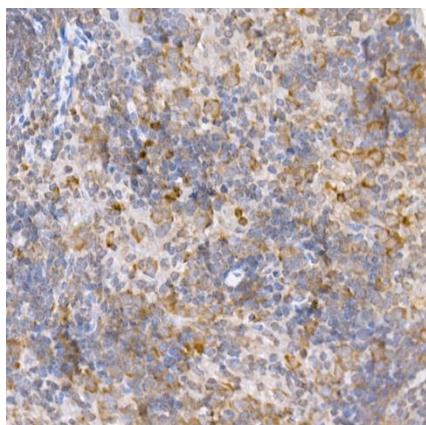
Immunohistochemistry of paraffin-embedded rat spleen using TRAF2 Polyclonal Antibody at dilution of 1:250 (40x lens).
 Performed high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded mouse brain using TRAF2 Polyclonal Antibody at dilution of 1:250 (40x lens).
 Performed high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded rat testis using TRAF2 Polyclonal Antibody at dilution of 1:250 (40x lens).
 Performed high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.



Immunohistochemistry of paraffin-embedded mouse spleen using TRAF2 Polyclonal Antibody at dilution of 1:250 (40x lens).
 Performed high-pressure antigen retrieval with 10 mM citrate buffer pH 6.0 before commencing with IHC staining protocol.