

Anti-TRAF2

Rabbit Polyclonal Antibody

Catalog # **T57-363BR**

Lot # O2109-46

Cited Applications

WB, ELISA, IF, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the TRAF2 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

TRAF2 antibody was raised against a 17 amino acid synthetic peptide from near the carboxy terminus of human TRAF2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

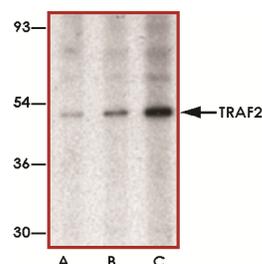
Scientific Background

TRAF2 (also known as TNF receptor-associated factor 2) belongs to the TNF receptor associated factor (TRAF) protein family that mediates signal transduction from members of the TNF receptor superfamily (1). TRAF2 is involved in TNF-alpha-mediated activation of MAPK8/JNK and NF-κB pathway. TRAF2 interacts with TNF receptors and forms a heterodimeric complex with TRAF1. The TRAF1-TRAF2 complex interacts with IAPs to mediate the anti-apoptotic signals from TNF receptors. TRAF2 also interacts with TRADD, a TNF receptor associated apoptotic signal transducer, to ensure the recruitment of IAPs for the direct inhibition of caspase activation (2).

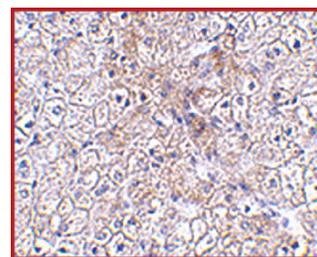
References

1. Wajant, H. Et al: The TNF-receptor-associated factor family: scaffold molecules for cytokine receptors, kinases and their regulators.". *Cell. Signal.* 2001;13 (6): 389-400.
2. Bradley, J.R., Poher, J.S. Tumor necrosis factor receptor-associated factors (TRAFs)". *Oncogene*,2001; 20 (44): 6482-9.

Sample Data



Western blot analysis of TRAF2 in human liver tissue lysate with TRAF2 antibody at (A) 0.5, (B) 1 and (C) 2 µg/ml.



Immunohistochemistry of TRAF2 in human liver tissue with TRAF2 antibody at 2.5 µg/ml.

Anti-TRAF2

Rabbit Polyclonal Antibody

Catalog Number

T57-363BR

Specific Lot Number

O2109-46

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.