

TRAF5 antibody (mAb)

Catalog No: 40920

RRID: AB_2793439

Clone: 55A219

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Concentration: 0.5 µg/µl

Molecular Weight: 66 kDa

Background: TRAF5 (TNF receptor-associated factor 5, RNF84) scaffold protein is a member of the tumor necrosis factor receptor-associated factor (TRAF) protein family and contains a meprin and TRAF homology (MATH) domain, a RING-type zinc finger, and two TRAF-type zinc fingers. TRAF proteins are associated with, and mediate signal transduction from, members of the TNF receptor superfamily. TRAF5 is one of the components of a multiple protein complex that binds to tumor necrosis factor (TNF) receptor cytoplasmic domains and mediates TNF-induced activation. TRAF5 mediates activation of NFκB (NFκB p50 & NFκB p65) and probably JNK. It also appears to be involved in apoptosis.

Immunogen: This TRAF5 antibody was raised against a fusion protein corresponding to amino acid residues 77 to 186 of human TRAF5.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

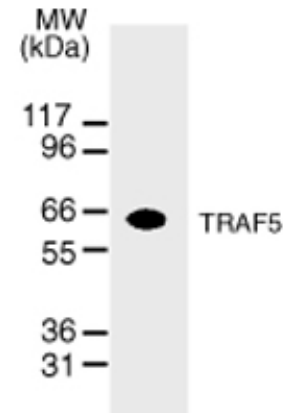
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TRAF5 mAb tested by Western blot.

TRAF5 detection by Western blot. The analysis was performed using 10 µg HeLa nuclear extract and TRAF5 mAb at a 2 µg/ml dilution.