

MALT1 antibody (mAb)

Catalog Nos: 39395, 39396

RRID: AB_2793240

Isotype: IgG

Application(s): WB

Reactivity: Human

Quantities: 200 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 92 kDa

Background: MALT1 enhances Bcl10-induced activation of NFκB and is also involved in nuclear export of Bcl10. MALT1 binds to TRAF6, inducing TRAF6 oligomerization and activation of its ligase activity. MALT1-dependent Bcl10 cleavage plays an important role in T-cell antigen receptor-induced integrin adhesion. MALT1 was independently identified as a protein of 824 amino acids, and human MALT2 isoform b is an amino acid protein of 813 amino acids. MALT1 is thought to play a role in NFκB signaling by enhancing NFκB (NFκB p50 & NFκB p65) activation through interaction with Bcl10. Interaction between MALT1 and Bcl10 mediates IKK (IKKα, IKKβ, IKKi/IKKε and IKKγ) activation *in vitro* through facilitating the ubiquitination of NEMO by the ubiquitin conjugating enzyme UBC13.

Immunogen: This MALT1 antibody was raised against a recombinant protein corresponding to amino acids 701-808 of mouse MALT1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

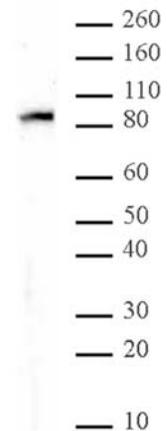
Application Notes:

Applications Validated by Active Motif:

WB: 0.5 - 2 µg/ml dilution

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.



MALT1 mAb tested by Western blot.

MALT1 detection by Western blot. The analysis was performed using 20 µg Jurkat nuclear extract and MALT1 mAb at a 2 µg/ml dilution.