

Caspase-14 antibody (mAb)

Catalog No: 40932

RRID: AB_2793447

Clone: 70A1426

Application(s): WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 30 kDa

Background: Caspase-14 (CASP14 or apoptosis-related cysteine peptidase 14) is thought to be a non-apoptotic caspase involved in epidermal differentiation. Caspase-14 has been shown to be processed and activated by Caspase-8 and Caspase 10 *in vitro*, and also by TNF-Related Apoptosis-Inducing Ligand (TRAIL) *in vivo*.

Immunogen: This Caspase-14 antibody was raised against a synthetic peptide corresponding to amino acid residues 2-18 of mouse Caspase-14.

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

Application Notes:

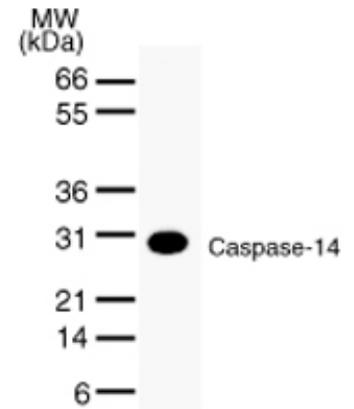
Applications Validated by Active Motif:

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. Store at 4°C for short term. 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.



Caspase-14 mAb tested by Western blot. Caspase-14 detection by Western blot. The analysis was performed using 10 µg of mouse spleen cell extract and Caspase-14 mAb at a 2 µg/ml dilution.