

Caspase-8 antibody (mAb)

Catalog No: 40930

RRID: AB_2793445

Clone: 90A992

Isotype: IgG1

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Concentration: 0.5 µg/µl

Molecular Weight: 55 kDa

Background: Caspase-8 (CASP8 or apoptosis-related cysteine peptidase 8) is the most upstream protease of the activation cascade of caspases responsible for the TNFRSF6/FAS mediated and TNFRSF1A induced cell death. Binding to the adaptor molecule FADD recruits it to either receptor. Caspase-8 cleaves and activates Caspase-3, Caspase-4, Caspase-6, Caspase-7, Caspase-9 and Caspase-10.

Immunogen: This Caspase-8 antibody was raised against a synthetic peptide corresponding to amino acid residues 217-222 of human Caspase-8.

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

Application Notes:

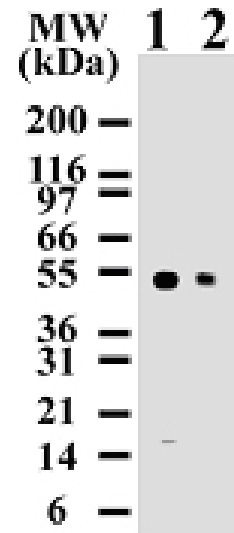
Applications Validated by Active Motif:

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.



Caspase-8 mAb tested by Western blot.

Caspase-8 detection by Western blot. The analysis was performed using Jurkat extracts with Caspase-8 mAb at 2 µg/ml (lane 1) and 0.5 µg/ml (lane 2) dilutions. The antibody only detects 55 kDa caspase-8 in staurosporine-treated (2 hours) Jurkat cells.