

Caspase-9 antibody (pAb) (Pro-form)

Catalog No: 40931

RRID: AB_2793446

Isotype: IgG

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.5 µg/µl

Molecular Weight: 46 kDa

Background: Caspase-9 (CASP9 or apoptosis-related cysteine peptidase 9) (Pro-form) is involved in the activation cascade of caspases responsible for apoptosis execution. Caspase-9 is processed by caspase APAF1, which is thought to be one of the earliest events in the caspase activation cascade. Caspase-9 is then involved in processing of Caspase-3.

Immunogen: This Caspase-9 antibody (Pro-form) was raised against a peptide corresponding to amino acid residues 299 -318 of human Pro-Caspase-9.

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

Application Notes:

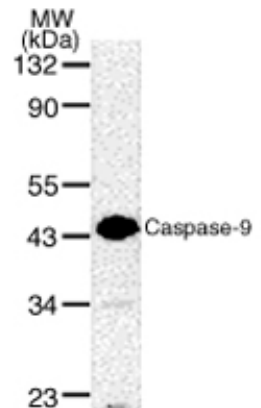
Applications Validated by Active Motif:

WB: 1:1,000 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-9 pAb (Pro-form) tested by Western blot.

Caspase-9 detection by Western blot. The analysis was performed using 50 µg HeLa whole-cell extract and Caspase-9 pAb (Pro-form) at a 1:1,000 dilution.