

## Caspase-7 antibody (mAb)

**Catalog No:** 40929

**RRID:** AB\_2793444

**Clone:** 25B204.1

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 35 kDa

**Background:** Caspase-7 (CASP7 or apoptosis-related cysteine peptidase 7) is involved in the activation cascade of caspases responsible for apoptosis execution. The precursor form of Caspase-7 is cleaved by Caspase-3 and Caspase-10, which is activated by cell death stimuli and induces apoptosis. Caspase-7 cleaves and activates sterol regulatory element binding proteins (SREBPs) and proteolytically cleaves poly(ADP-ribose) polymerase (PARP, see PARP-1 C-terminal and PARP-1 N-terminal antibodies).

**Immunogen:** This Caspase-7 antibody was raised against a recombinant human Caspase-7 protein.

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

### Application Notes:

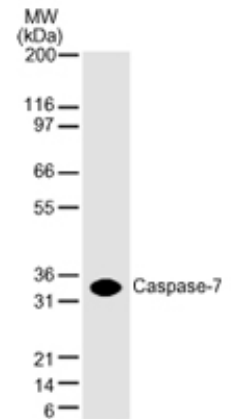
Applications Validated by Active Motif:

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-7 mAb tested by Western blot.

Caspase-7 detection by Western blot. The analysis was performed using Jurkat nuclear extracts and Caspase-7 mAb at dilutions of 2 µg/ml (lane 1) and 0.5 µg/ml (lane 2). This antibody only detects the 35 kDa caspase-7 corresponding to pro-caspase-7 in non-apoptotic Jurkat cells.