

## A20 antibody (mAb)

**Catalog No:** 40901

RRID: AB\_2793426

Clone: 59A426

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 1 µg/µl

Molecular Weight: 70 kDa

**Background:** A20 is a zinc finger protein induced by Tumor Necrosis Factor (TNF) and involved in limiting inflammation through inhibition of NFκB (NFκB p50 & NFκB p65) activation. A20 also inhibits TNF-mediated apoptosis. A20 is also known as TNFA1P2, TNFAIP3 and OTUD7C.

**Immunogen:** This A20 antibody was raised against a full-length protein corresponding to human A20.

**Buffer:** PBS containing 0.05% BSA and 0.05% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

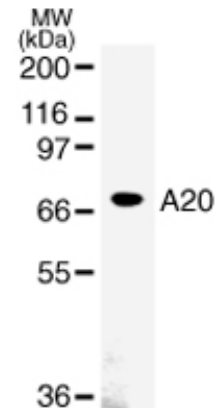
Applications Validated by Active Motif:

WB: 3 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.

**A20 mAb tested by Western blot.**

Detection of A20 by Western blot. A20 can be detected in Daudi whole-cell extract at 3 µg/ml.