

DAP-3 antibody (mAb)

Catalog No: 40963

RRID: AB_2793460

Clone: 42C617

Application(s): WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 45 kDa

Background: DAP-3 (Death Associated Protein 3) potentially binds ATP/GTP and may be a functional partner of the mitoribosomal protein S27. The DAP-3 gene encodes a 28S subunit protein that also participates in apoptotic pathways that are initiated by tumor necrosis factor-alpha, Fas ligand, and gamma interferon.

MW
116
97

Immunogen: This DAP-3 antibody was raised against a synthetic peptide corresponding to amino acid residues 51-66 of human DAP-3.

66
55

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

36
31
21
14

 **DAP-3**

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.

DAP-3 mAb tested by Western blot.

DAP-3 detection by Western blot. The analysis was performed using 10 µg Jurkat whole-cell extract and the DAP-3 mAb at a 1 µg/ml dilution.

This product is for research use only and is not for use in diagnostic procedures.