

DAP-5 antibody (mAb)

Catalog No: 40964**RRID:** AB_2793461**Clone:** 39C534**Application(s):** WB**Reactivity:** Human, Mouse**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Mouse**Isotype:** IgG1**Concentration:** 0.5 µg/µl**Molecular Weight:** 97 kDa

Background: DAP-5 (eIF4G2 or eukaryotic translation initiation factor 4 gamma 2) shares similarity with the C-terminal region of eIF4G, which contains the binding sites for eIF4A and eIF3. eIF4G, in addition, contains a binding site for eIF4E at the N-terminus. Unlike eIF4G, which supports cap-dependent and independent translation, this gene product functions as a general repressor of translation by forming translationally inactive complexes. *In vitro* and *in vivo* studies indicate that translation of this mRNA initiates exclusively at a non-AUG (GUG) codon. Alternatively spliced transcript variants encoding different isoforms of this gene have been described. DAP-5 appears to play a role in the switch from cap-dependent to IRES-mediated translation during mitosis, apoptosis and viral infection.

Immunogen: This DAP-5 antibody was raised against a synthetic peptide corresponding to amino acid residues 796-812 of human DAP-5.

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

Application Notes:

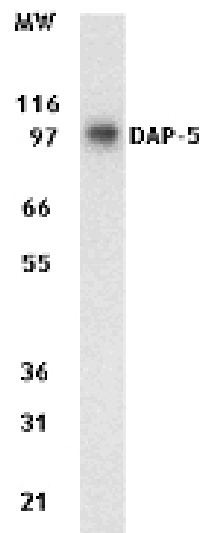
Applications Validated by Active Motif:

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. 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.

**DAP-5 mAb tested by Western blot.**

DAP-5 detection by Western blot. The analysis was performed using 15 µg of Jurkat whole-cell extract and the DAP-5 mAb at a 1 µg/ml dilution.