

DRAK1 antibody (pAb)

Catalog No: 40923

RRID: AB_2793440

Application(s): WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Concentration: 0.5 µg/µl

Molecular Weight: 56 kDa

Background: Overexpression of DRAK1 (STK17A or DAP kinase-related apoptosis-inducing protein kinase 1) and ZIP kinase can cause the morphological changes typical of apoptosis in NIH/3T3 cells. Both DRAKs and ZIP kinase share significant homology at the amino acid level. The kinase domains of both DRAKs and ZIP kinase are homologous to DAP (Death-associated protein) kinase, which is involved in the apoptotic signaling induced by interferon-γ. Overexpression of DRAKs in NIH/3T3 cells lead to apoptosis. These experiments also indicate that the C-terminal domain of DRAKs are important for induction of apoptosis.

Immunogen: This DRAK1 antibody was raised against a peptide corresponding to amino acid residues 5 to 9 of human DRAK1.

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

Application Notes:

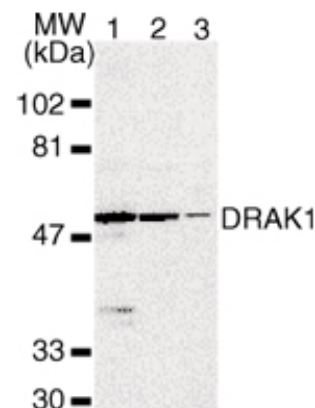
Applications Validated by Active Motif:

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.



DRAK1 pAb tested by Western blot.

Detection of DRAK1 by Western blot. The analysis was performed using MOLT4 (lane 1), A-431 (lane 2) and NIH/3T3 (lane 3) nuclear extracts and DRAK1 pAb at a 2 µg/ml dilution.