

DR6 antibody (pAb)

Catalog No: 40927

RRID: AB_2793443

Isotype: IgG

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 66 kDa

Background: DR6 (also known as Tumor Necrosis Factor Receptor Superfamily, member 21, TNFRSF21) is a member of the TNF-receptor superfamily. This receptor interacts with TRADD to mediate signal transduction of TNF receptors. It may activate NFκB (NFκB p50 & NFκB p65) and MAPK8/JNK and induce apoptosis.

Immunogen: This DR6 antibody was raised against a cytoplasmic domain corresponding to amino acid residues 371-655 of human DR6.

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

Application Notes:

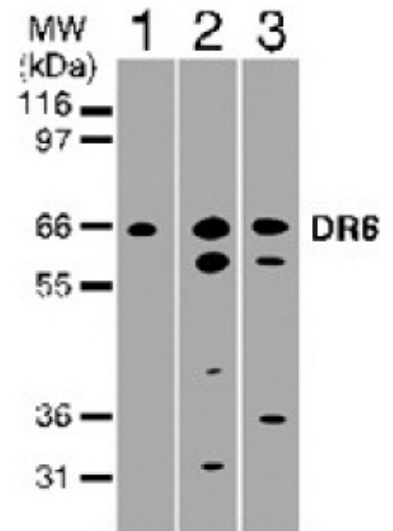
Applications Validated by Active Motif:

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.



DR6 pAb tested by Western blot analysis.

Western blot analysis of 30 µg of total cell lysate from Jurkat (lane 1), HeLa (lane 2), and Daudi (lane 3) cells tested at 2 µg/ml.