

DR4 antibody (mAb)

Catalog No: 40934

RRID: AB_2793449

Clone: 32A1380

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Isotype: IgG1

Concentration: 0.5 µg/µl

Molecular Weight: 57 kDa

Background: DR4 (also known as Tumor Necrosis Factor Receptor Superfamily, member 10a, TNFRSF10A, or TRAILR1) is a member of the TNF-receptor superfamily that is activated by TRAIL and induces apoptosis.

Immunogen: This DR4 antibody was raised against a peptide corresponding to amino acid residues 1-20 of human DR4.

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

Application Notes:

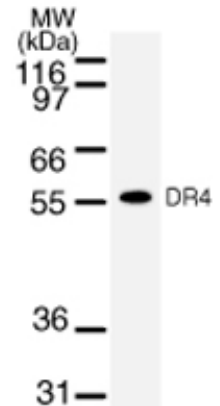
Applications Validated by Active Motif:

WB: 2 - 4 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.



DR4 mAb tested by Western blot.

Detection of DR4 by Western blot. The analysis was performed using Daudi whole-cell extract and DR4 mAb at a 2 µg/ml dilution.