

## DcR3 antibody (mAb)

**Catalog No:** 40933

**RRID:** AB\_2793448

**Clone:** 37A565

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 34 kDa

**Background:** DcR3 (TNFRSF6B or tumor necrosis factor receptor superfamily member 6b) is thought to play a regulatory role in suppressing FasL- and LIGHT-mediated cell death. DcR3 acts as a decoy receptor that competes with death receptors for ligand binding. Overexpression of this gene has been noted in gastrointestinal tract tumors, and it is located in a gene-rich cluster on chromosome 20, with other potentially tumor-related genes. Two transcript variants encoding the same isoform, but differing in the 5' UTR, have been observed for this gene.

**Immunogen:** This DcR3 antibody was raised against a full-length protein corresponding to human DcR3.

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

### Application Notes:

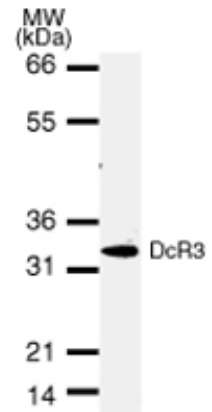
Applications Validated by Active Motif:

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.



### DcR3 mAb tested by Western blot.

Detection of DcR3 by Western blot. The analysis was performed using 30 µg Jurkat whole-cell extract and DcR3 mAb at a 2 µg/ml dilution.