

## RANK antibody (mAb)

**Catalog No:** 40917

**RRID:** AB\_2793437

**Clone:** 9A725

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Mouse

**Isotype:** IgG1

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 97 kDa

**Background:** RANK (TNFRSF11A, CD265, PDB2, TRANCER) can interact with various TRAF family proteins (TRAF2, TRAF5), through which this receptor induces the activation of NFκB and MAPK8/JNK. RANK and its ligand are important regulators of the interaction between T cells and dendritic cells. RANK is also an essential mediator for osteoclast and lymph node development. Studies of the mouse counterpart suggest that this receptor directly mediates the osteoprotegerin (OPG) ligand (OPGL)-induced osteoclastogenesis in osteoclast precursor cells.

**Immunogen:** This RANK antibody was raised against a GST-fusion protein containing amino acid residues 326-616 of human RANK.

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

### Application Notes:

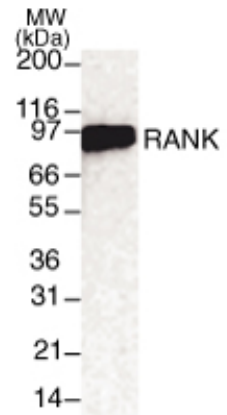
Applications Validated by Active Motif:

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.



### RANK mAb tested by Western blot.

RANK detection by Western blot. The analysis was performed using 293 cells transfected with RANK cDNA and RANK mAb at a 2 µg/ml dilution. A single protein band of approximate mol. wt. of 97 kDa was detected.