

## TRANCE/RANKL antibody (mAb)

**Catalog No:** 40940

**Clone:** 12A380

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human, Mouse

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 0.5 µg/µl

**Quantity:** 100 µg

**Molecular Weight:** 40 kDa

**Background:** TRANCE / RANKL – Tumor Necrosis Factor-Related Activation-induced Cytokine, also known as Receptor Activator of NFκB Ligand is a member of the tumor necrosis factor (TNF) cytokine family and is a ligand for osteoprotegerin (OPG). TRANCE/RANKL functions as a key factor for osteoclast differentiation and activation, is a key survival factor for dendritic cells and is involved in the regulation of T cell-dependent immune response. Overexpression of this protein leads to an increase of osteoclastogenesis and bone loss. This protein may also have a role in regulating apoptosis, by activating the anti-apoptotic kinase AKT/PKB through a signaling complex involving SRC kinase and tumor necrosis factor receptor-associated factor 6 (TRAF6).

**Immunogen:** This TRANCE/RANKL antibody was raised against a GST-fusion protein containing amino acid residues 1-317 of human TRANCE/RANKL.

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

**Application Notes:**

Validated Applications:

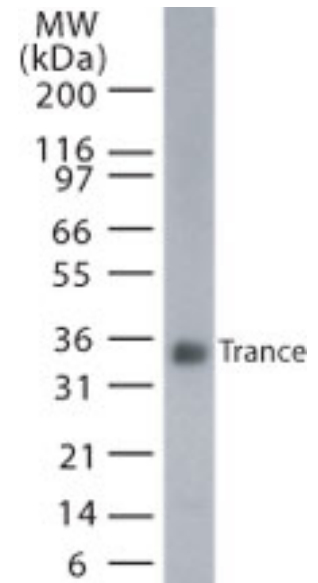
WB: 4 - 6 µg/ml dilution

Published Applications: see references

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all Blotto 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. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**TRANCE/RANKL mAb tested by Western blot.**

TRANCE detection by Western blot. The analysis was performed using TRANCE/RANKL mAb and transfected RAW cells.