## RANKL monoclonal antibody (12A380)

RANKL (TRANCE, OPGL, ODF) is a member of the TNF family. RANKL (receptor activator of NF-kB ligand) or TRANCE (TNF-related activation-induced cytokine receptor) has been implicated in interactions between T cells and dendritic cells. RANKL binds to RANK on dendritic cells, upregulates the expression of anti-apoptotic protein BcL-X<sub>L</sub> suggesting a role in dendritic cell survival. RANKL is also important in T- and B-cell maturation. As OPGL (osteoprotegerin ligand) or ODF (osteoclast differentiation factor), the same protein can both activate mature osteoclasts and mediate osteoclastogenesis. RANKL deficient mice show severe osteoporesis and complete abscence of osteoclasts as a result of lack of osteogenesis.

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

**Ordering Information** 

**Order Online** »

ALX-804-244-C100

100µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Use/Stability** Stable for at least 6 months when stored at +4°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name ODF, TRANCE, OPGL, TNFSF 11, CD254

**Application** Flow Cytometry, IHC (FS), WB

**Application Notes**Detects a band of ~40kDa by Western blot.

**Clone** 12A380

**Formulation** Liquid. In PBS containing 0.05% BSA and 0.05% sodium

azide.

**Host** Mouse

**Immunogen** Recombinant mouse RANKL (aa 1-317).

lsotype IgG1

Positive Control Activated EL4 cells (mouse) or Saos-2 (human) cells can

be used as positive control.

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Western Blot (4-6µg/ml)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID O35235

Last modified: May 29, 2024



info-

eu@enzolifesciences.com