

RANKL monoclonal antibody (12A380)

RANKL (TRANCE, OPGL, ODF) is a member of the TNF family. RANKL (receptor activator of NF- κ B 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_L 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 absence 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

100 μ g

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).
Isotype	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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.