

# RANKL monoclonal antibody (12A380)

RANKL (TRANCE, OPGL, ODF) is a member of the TNF family. RANKL (receptor activator of NF- $\kappa$ 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<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 absence of osteoclasts as a result of lack of osteogenesis.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months when stored at +4°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	ODF, TRANCE, OPGL, TNFSF 11, CD254
<b>Application</b>	Flow Cytometry, IHC (FS), WB
<b>Application Notes</b>	Detects a band of ~40kDa by Western blot.
<b>Clone</b>	12A380
<b>Formulation</b>	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant mouse RANKL (aa 1-317).
<b>Isotype</b>	IgG1
<b>Positive Control</b>	Activated EL4 cells (mouse) or Saos-2 (human) cells can be used as positive control.
<b>Purity Detail</b>	Protein G-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (4-6µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human, Mouse
<b>UniProt ID</b>	O35235

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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