

RANKL monoclonal antibody (12A668)

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-2430	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

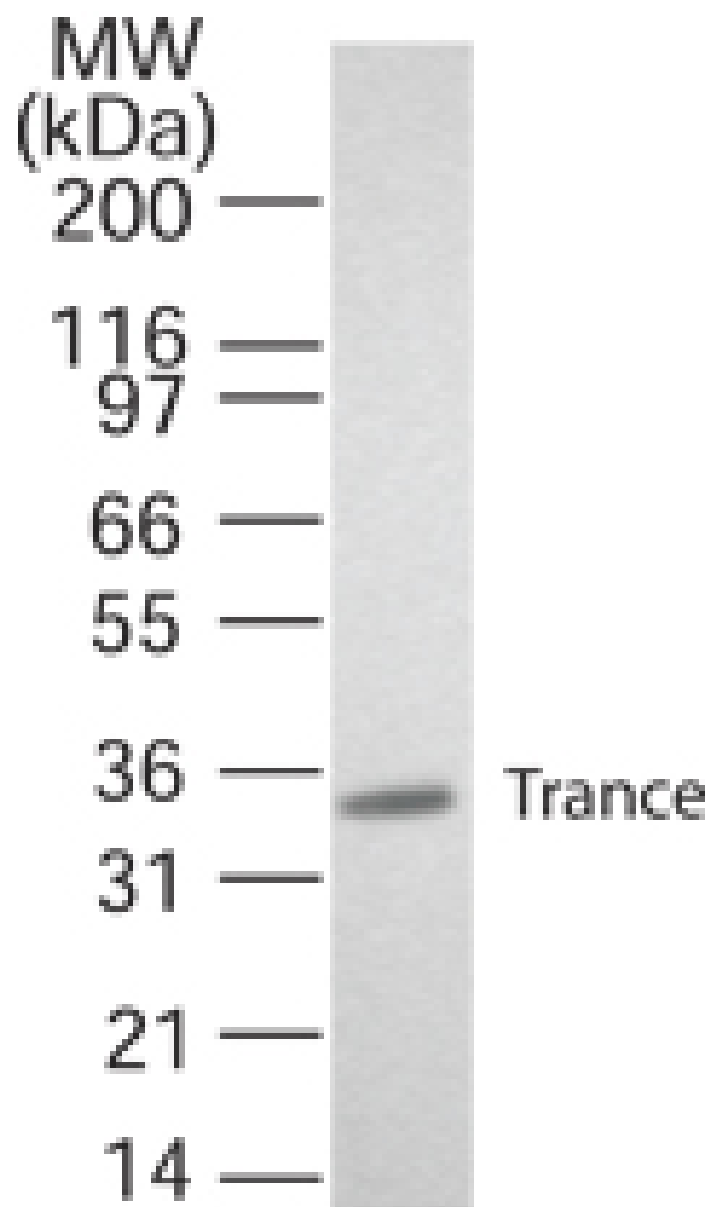
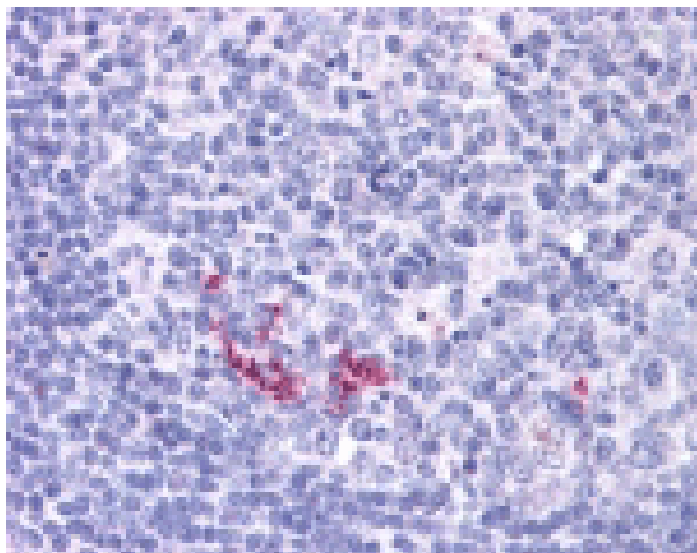


Figure 1: Western blot analysis of RANKL in transfected cells using MAb to RANKL (12A668) (Prod. No. ALX-804-243).

Handling & Storage

Use/Stability	Stable for 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	IHC (PS), WB
Application Notes	Detects a band of ~35-40kDa by Western blot. Not suitable for flow cytometry.
Clone	12A668
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant mouse RANKL (aa 1-317).
Isotype	IgG1 κ
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 5 μ g/ml)Western Blot (1-2 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
Technical Info / Product Notes	Replacement for ADI-AAM-425 , see also ADI-AAM-425AF

UniProt ID

O35235

Worry-free Guarantee

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

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)