## RANKL monoclonal antibody (12A668)

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

**Ordering Information** 

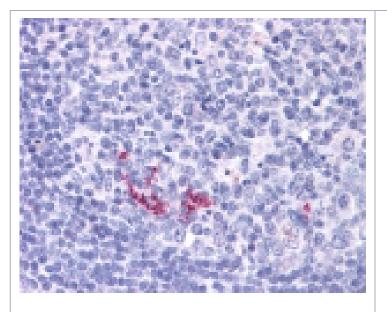
Order Online »

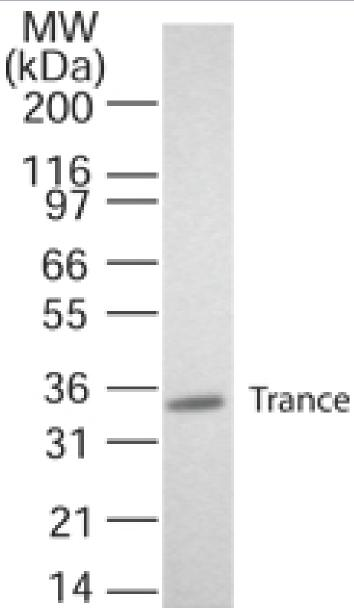
ALX-804-243-C100

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 ĸ

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 5µg/ml)Western

Blot (1-2µg/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

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



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

eu@enzolifesciences.com

Phone: 800.942.0430

info-

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

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

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

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