

RANKL (human) monoclonal antibody (Ranky-1)

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

Ordering Information

[Order Online »](#)

ALX-804-830-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

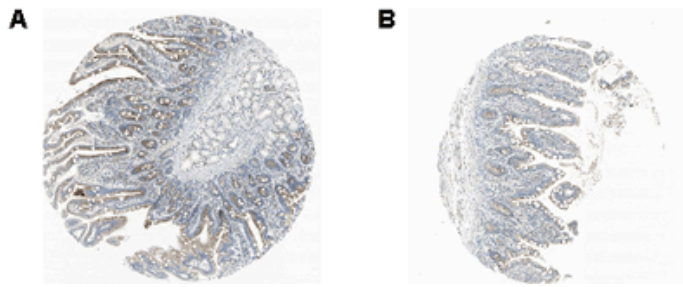


Figure: Immunohistochemical staining of endogenous human RANKL in different human tissues (paraffin sections) using MAb to RANKL (human)(Ranky-1) (Prod. No. [ALX-804-830](#)). **Method:** Different human tissues (A: Duodenum; B: Small intestine) were stained with MAb to RANKL (human) (Ranky-1) by standard immunohistochemistry. *Pictures courtesy of Human Protein Atlas ([www.proteinatlas.org](#))*. ⁶⁶⁷⁷⁷⁷⁷⁷

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
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	ELISA, IHC (PS), WB
Application Notes	Does not work for Flow Cytometry or Immunoprecipitation. Does not block RANK-RANKL interaction.
Clone	Ranky-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human soluble RANKL extracellular domain (aa 151-316) (Prod. No. ALX-522-012).
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (capture)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human

UniProt ID

O14788

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)