

# IL-33 monoclonal antibody (Nessy-1) (preservative free)

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

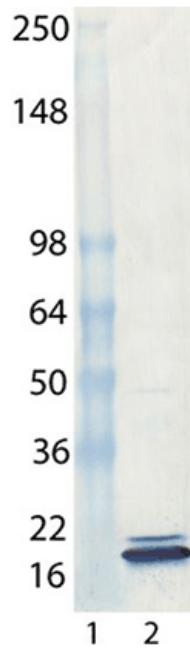
## Ordering Information

[Order Online »](#)

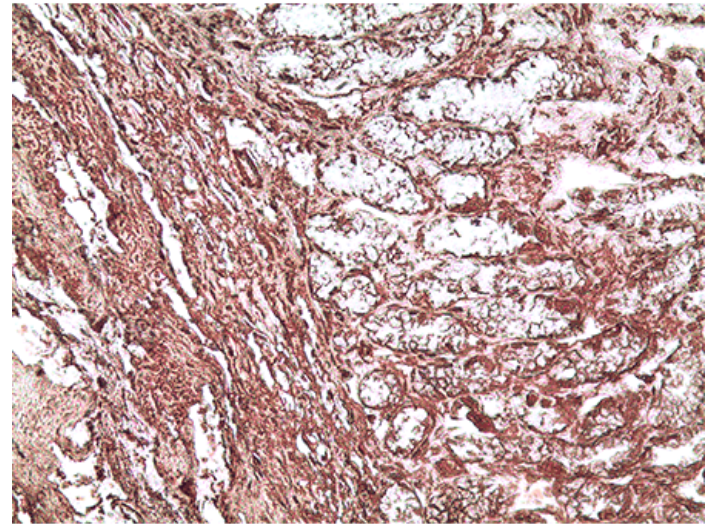
ENZ-ABS665-0100	Preservative Free - 100µg
-----------------	---------------------------

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS665: Lane 1: MW Marker, Lane 2: IL-33 (human), (recombinant) (Prod. No. ALX-522-098).



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human stomach tissue stained with IL-33 monoclonal antibody (Nessy-1), (Prod. No. ENZ-ABS665) at 10µg/ml.

## 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	Interleukin-33, IL-1F11, NF-HEV
Application	ELISA, IHC (PS), IP, WB
Clone	Nessy-1
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Recombinant human IL-33 (aa 112-270) (Prod. No. ALX-522-098).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500 – 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 hybridoma tissue culture supernatant.
Species Reactivity	Human, Mouse, Rat
UniProt ID	O95760

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)