

IL-33 monoclonal antibody (Nessy-1) (biotin conjugate)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-840B-C100	Biotin - 100µg
-------------------	----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
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)
Clone	Nessy-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Gene/Protein Identifier	90865 (Entrez GeneID), 608678 (OMIM)
Host	Mouse
Immunogen	Recombinant human IL-33 (aa 112-270) (Prod. No. ALX-522-098).
Isotype	IgG1κ
Purity Detail	Protein G-affinity purified.
Source	Purified from ascites.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes pro-IL-33 (precursor) and mature IL-33.
UniProt ID	O95760
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



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)