

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

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|--------------------|---------------------------|
| ALX-804-840PF-C100 | Preservative Free - 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), IP, WB |
| Clone | Nessy-1 |
| Formulation | Liquid. In PBS. |
| 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. |
| Recommendation Dilutions/Conditions | Immunoprecipitation (1:100)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 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)