

IL-33 polyclonal antibody (AT110)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-933-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

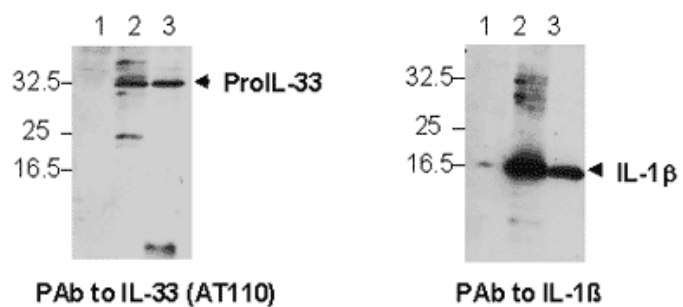


Figure: Endogenous IL-33 is detected in the supernatant of THP1 using PAb to IL-33 (AT110) (Prod. No. [ALX-210-933](#)). **Method:** THP-1 cells were treated with PBS (lane 1), monosodium urate (MSU) (Prod. No. ALX-400-047) (lane 2) or peptidoglycan (PGN) (lane 3). The supernatants were collected

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store 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, NF-HEV, IL-1F11
Application	ELISA, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Gene/Protein Identifier	90865 (Entrez GeneID), 608678 (OMIM)
Host	Rabbit
Immunogen	Recombinant human soluble IL-33 (interleukin-33) (aa 112-270) (Prod. No. ALX-522-098).
Positive Control	Prod. No. ALX-522-098; Prod. No. ALX-522-101
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	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 rabbit serum.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes pro-IL-33 and mature IL-33.

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)