

# IL-33 polyclonal antibody (AT110)

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

Citations: 4

[View Online »](#)

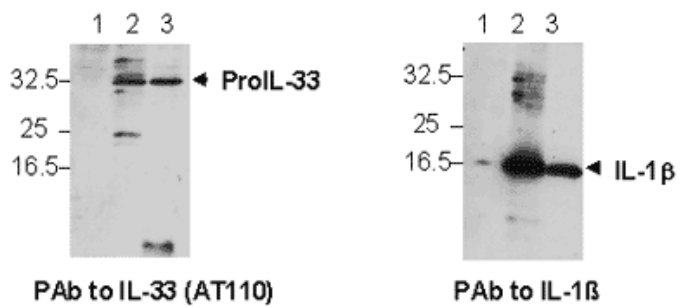
Ordering Information

[Order Online »](#)

ALX-210-9330	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

<b>Use/Stability</b>	Stable for at least 1 year after receipt when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Interleukin-33, NF-HEV, IL-1F11
<b>Application</b>	ELISA, WB
<b>Formulation</b>	Liquid. In PBS containing 0.02% sodium azide.
<b>Gene/Protein Identifier</b>	90865 (Entrez GeneID), 608678 (OMIM)
<b>Host</b>	Rabbit
<b>Immunogen</b>	Recombinant human soluble IL-33 (interleukin-33) (aa 112-270) (Prod. No. ALX-522-098).
<b>Positive Control</b>	Prod. No. ALX-522-098; Prod. No. ALX-522-101
<b>Purity Detail</b>	Protein A-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from rabbit serum.
<b>Species Reactivity</b>	Human, Mouse, Rat
<b>Specificity</b>	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)