

IL-33 (mouse), (recombinant)

IL-33 is a member of the IL-1 family. It is involved in T helper type 2 response via the activation of its receptor ST2 (IL-1RL1) driving the production of Th2-associated cytokines. The stimulation of ST2 by IL-33 induces the NF-κB and MAPK pathways. The mouse IL-33 protein is expressed with a prodomain (aa 1-108) which is cleaved to produce the 18kDa mature protein (109-266). Historical data has shown that IL-33 activates mouse and human ST2-dependent NF-κB pathway and induces IL-6 production in P815 mastocytoma cells and in bone marrow derived mast cells.

Citations: 7

[View Online »](#)

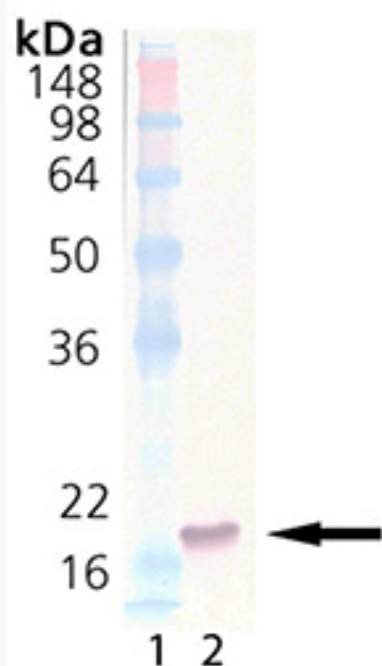
Ordering Information

[Order Online »](#)

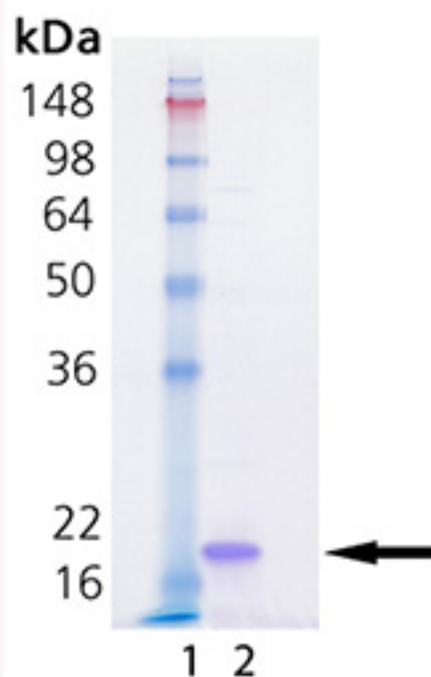
ALX-522-101-C010	10µg
ALX-522-101-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis: Lane 1: MW Marker; Lane 2: 100ng of purified IL-33 (mouse), (recombinant) protein probed with IL-33 mAb (Prod. no. ALX-804-840PF-C100) at 0.5ug/ml.



SDS-PAGE analysis: Lane 1: MW Marker; Lane 2: 1.0ug of purified IL-33 (mouse), (recombinant).

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°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
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~18kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile distilled water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in <i>E. coli</i> . Untagged mouse IL-33 (aa 109-266).
Specificity	Binds to mouse and human ST2.
UniProt ID	Q8BVZ5

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)