

IL-33 monoclonal antibody (Nessy-1)

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

Citations: 63

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-840-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

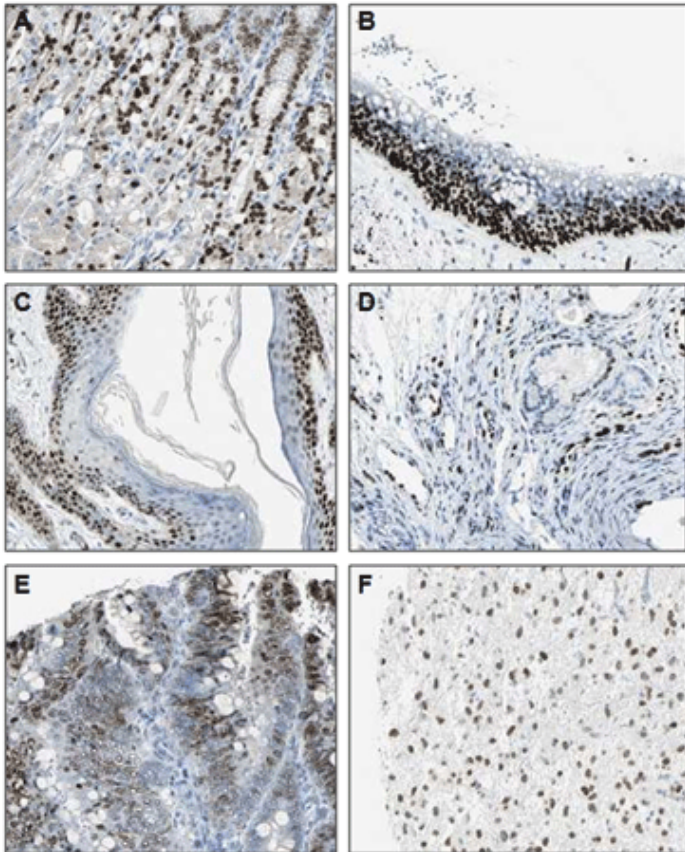
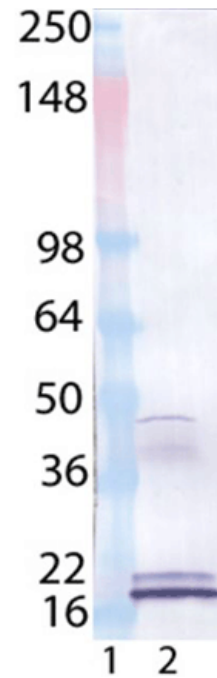
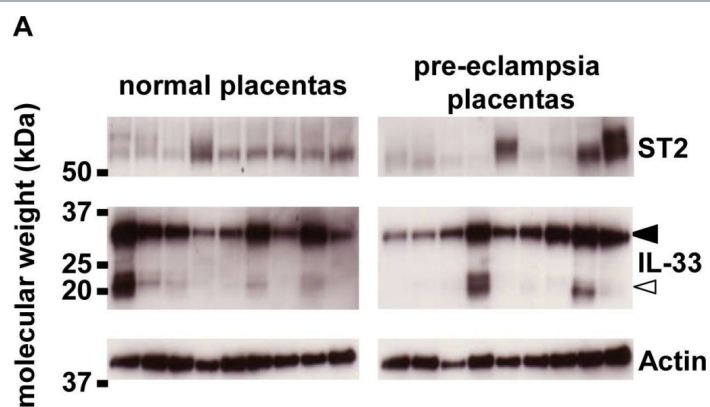


Figure: Immunohistochemical staining of endogenous human IL-33 in different human tissues (paraffin sections) using MAb to IL-33 (Nessy-1) (Prod. No. ALX-804-840). Method: Different human normal tissues (A: Stomach; B: Bronchus; C: Vulva/Anal skin; D: Nasopharynx) or cancer tissues (E: Colorectal cancer; F: Malignant glioma) were stained with MAb to IL-33 (Nessy-1) by standard immunohistochemistry. Pictures courtesy of Human Protein Atlas (<http://www.proteinatlas.org/>). """"""

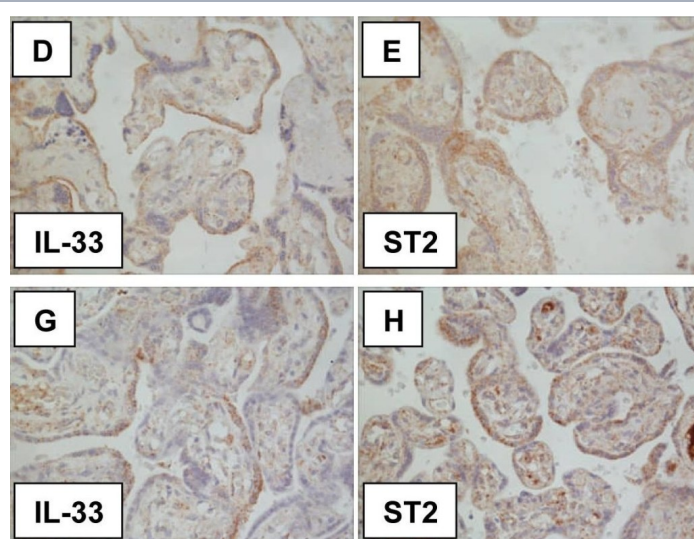


Western Blot analysis of ALX-804-840: Lane 1: MW Marker, Lane 2: IL-33 (human), (recombinant), (Prod. No. ALX-522-098).



Expression of IL-33 and sST2 by the placenta. A) IL-33 was detected by western blotting of placental lysates from normal and pre-eclampsia placentas (n = 9). Both the full length (closed arrow) and cleaved forms (open arrow) were detected. ST2 was detected in each of the placental lysates. B) Densitometry of protein bands standardised to actin expression revealed no change in the levels of IL-33 nor sST2 between normal and pre-eclamptic placentas. C–H) Expression of IL-33 and sST2 can be seen in by immunocytochemical analysis of both normal (n = 3) and pre-eclampsia (n = 3) placentas (C–E and F–H show staining from two representative placentas respectively).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: ST2 and IL-33 in pregnancy and pre-eclampsia. *PLoS One* (2011)



Expression of IL-33 and sST2 by the placenta. A) IL-33 was detected by western blotting of placental lysates from normal and pre-eclampsia placentas (n = 9). Both the full length (closed arrow) and cleaved forms (open arrow) were detected. ST2 was detected in each of the placental lysates. B) Densitometry of protein bands standardised to actin expression revealed no change in the levels of IL-33 nor sST2 between normal and pre-eclamptic placentas. C–H) Expression of IL-33 and sST2 can be seen in by immunocytochemical analysis of both normal (n = 3) and pre-eclampsia (n = 3) placentas (C–E and F–H show staining from two representative placentas respectively).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: ST2 and IL-33 in pregnancy and pre-eclampsia. *PLoS One* (2011)

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 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.
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.

Technical Info / Product Notes

Cited samples:

[For an overview on cited samples please click here.](#)

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)