

IL-33 monoclonal antibody (03)

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

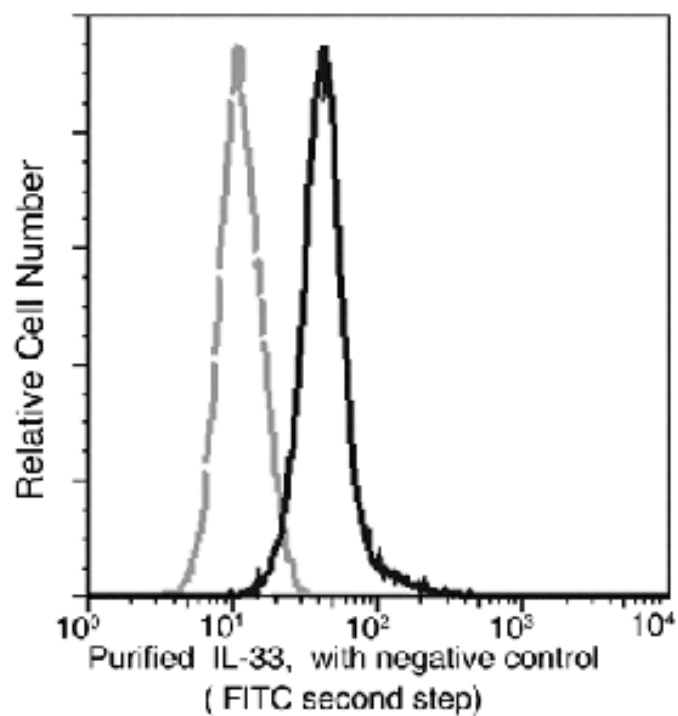
Ordering Information

[Order Online »](#)

ENZ-ABS948-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Purified anti-Human IL-33 on HUVEC cells. HUVEC cells were treated and stained with Purified Mouse anti-IL-33 (Bold line histogram), then stained with a FITC-conjugated second step antibody.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

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, Flow Cytometry

Clone 03

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Mouse

Immunogen Recombinant human IL-33 protein.

Isotype IgG2b

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000)Flow Cytometry (1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

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