

IL-33 (human), ELISA kit

The IL-33 (human), ELISA kit is a colorimetric immunometric immunoassay kit

The IL-33 (human), EIA kit is a colorimetric immunometric immunoassay kit with results in 3 hours.

IL-33, also named Nuclear Factor from High Endothelial Venules (NF-HEV), is a member of the Interleukin-1 (IL-1) family. Members of this family, such as IL-1 α / β and IL-18, are known to play important roles in host defense, immune regulation, neuronal injury and inflammation. IL-33 has been reported to be the specific ligand for ST2, an orphan receptor of the IL-1 receptor family. Stimulation of ST2 by IL-33 induces NF- κ B and MAPK signaling pathways. IL-33 is synthesized in response to cytokines such as IL-1 β and TNF- α and plays an important role in TH2-associated immunology by triggering increased secretion of IL-5 and IL-13 from polarized TH2 cells. IL-33 has also been found to be localized in the nucleus of human epithelial and endothelial cells, and can be present in serum and synovial fluid samples of patients with rheumatoid arthritis. Overexpression of IL-33 in mice has been found to induce asthmatic symptoms, and enhance joint inflammation.

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

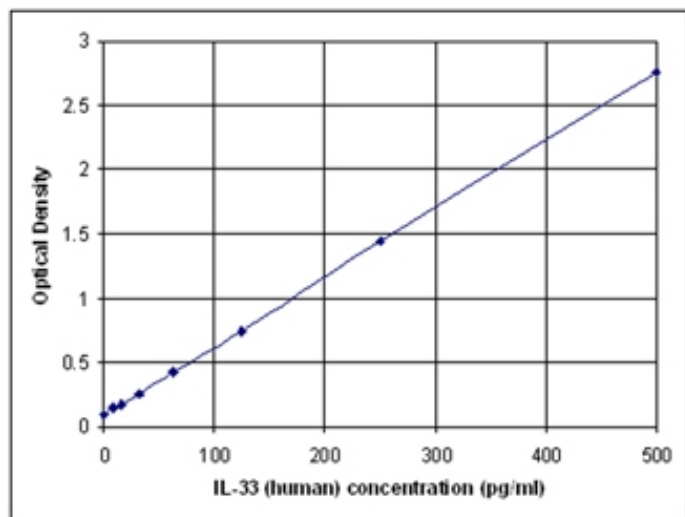
ADI-900-201

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - reliably measure as little as 1.7pg/ml of IL-33
- Low dynamic range provides sensitive measurements
- Specific - low cross-reactivity to other interleukins
- Non-radioactive - enzymatic assay format is safe and reliable
- Easy-to-use - liquid color-coded reagents reduce errors



Handling & Storage

Use/Stability	Store all components at +4°C, except Standard at -20°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin-33, NF-HEV
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of human IL-33 in cell lysates, culture supernatants, serum, plasma and synovial fluid samples.
Assay Time	3 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	Microtiter Plate, Conjugate, Antibody, Assay Buffer 13, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Extraction Reagent Concentrate
Crossreactivity	Human IL-33 (100%), mouse IL-33 (<0.02%). Unaffected by human IL-1 α , IL-2, IL-10, MCP-1, IL-13, MIP-1 β , IL-4, IL-6, IL-8, TNF α , IFN γ , IL-1 β IL-5, IL-15, IL-17, IL-21, IL-29, RANTES
Quantity	1 x 96-well plate
Sensitivity	1.7 pg/ml (range 7.8 – 500 pg/ml)
Species Reactivity	Human
Technical Info / Product Notes	Cited samples: Cytokine ELISAs Cited Samples

UniProt ID O95760

Wavelength 450 nm

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)