

IL-33 (human), (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 human IL-33 protein is expressed with a prodomain (aa 1-111) which is cleaved by caspase-1 to produce the 18kDa mature protein (112-270). Historical data has shown that IL-33 activates human ST2-dependent NF-κB pathway, but it does not activate mouse ST2-dependent NF-κB.

Citations: 3

[View Online »](#)

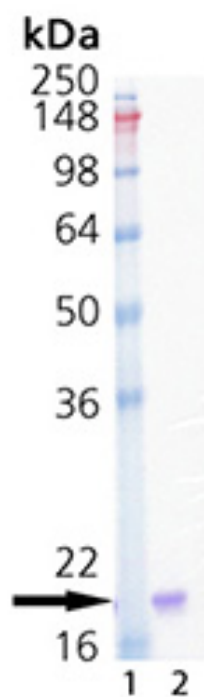
Ordering Information

[Order Online »](#)

ALX-522-098-C010	10µg
ALX-522-098-3010	SuperPack - 3x10µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MWM; Lane 2: 1.0 μg of purified IL-33 (human), (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 Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~18kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
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 human IL-33 (aa 112-270).
Specificity	Binds to human ST2.
UniProt ID	O95760