

IL-33 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat IL-33 (IN)

Instruction Manual

Catalog Number	PK-AB718-4273
Synonyms	IL-33 Antibody; Interleukin-33, Nuclear factor from high endothelial venules, NFHEV
Description	Interleukin-33 (IL-33) is a recently identified member of the IL-1 family of cytokines whose other members include IL-1a/b, IL-1Ra and IL-18. Its receptor has been shown to be ST2, an IL-1 receptor family member that also acts as a negative regulator of TLR-IL-1R signaling and IL-1R accessory protein (IL-1RAcP). Receptor binding of IL-33 activates NF-kappaB and MAP kinases and induces the expression of TH2-associated cytokines such as IL-4, IL-5 and IL-6. Prolonged IL-33 treatment of mice led to the development of eosinophilia, splenomegaly, and severe pathological changes in mucosal organs such as lungs, esophagus and small intestine. Recent experiments have shown that IL-33 can also co-localize with heterochromatin and possesses transcriptional repressor activities, indicating that IL-33 may function as both a proinflammatory cytokine and an intracellular nuclear factor with transcriptional regulatory properties. Despite its predicted molecular weight, IL-33 will often run at higher molecular weight in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	IL-33 antibody was raised in rabbits against a 19 amino acid peptide from near the center of human IL-33.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. IL-33 antibody can be used for the detection of IL-33 by Western blot at 1 - 2 µg/mL. Despite its predicted molecular weight, IL-33 will often run at higher molecular weight in SDS-PAGE. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Schmitz J, Owyang A, Oldham E, et al. IL-33, and interleukin-1-like cytokine that signals via the IL-1 receptor-related protein ST2 and induces T helper type 2-associated cytokines. <i>Immunity</i> 2005; 23:479-90. Dinarello CA. Interleukin-18, a proinflammatory cytokine. <i>Eur. Cytokine Netw.</i> 2000; 11:483-6. Brint EK, Xu D, Liu H, et al. ST2 is an inhibitor of interleukin 1 receptor and Toll-like receptor 4 signaling and maintains endotoxin tolerance. <i>Nat. Immunol.</i> 2004; 5:373-9. Chackerian AA, Oldham ER, Murphy EE, et al. IL-1 receptor accessory protein and ST2 comprise the IL-33 receptor complex. <i>J. Immunol.</i> 2007; 179:2551-5.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4273P; IL-33 Peptide Cat.No. PK-AB718-1369; Human Lymph Node Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.