

Product Information

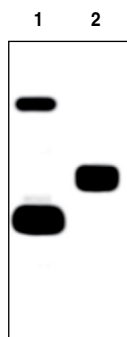


Interleukin-33 (human) Monoclonal Antibody (Clone IL33305B)

Item No. 10809

Contents:	This vial contains a 1 mg/ml solution of antibody in PBS, pH 7.4.
Synonyms:	IL-1F11, IL-33, NF-HEV
Antigen:	Recombinant human IL-33
Host:	Mouse, clone IL33305B
Isotype:	IgG _{2ak}
Cross Reactivity:	(+) Human IL-33; (-) mouse IL-33
Stability:	≥6 months at -20°C
Applications:	ELISA, immunohistochemistry (IHC), immunoprecipitation, and western blot (WB). The recommended starting dilutions are: ELISA (direct or indirect): 1:2,000-1:5,000, IHC: 1:50-1:200, and WB: 1:2,000-1:5,000.

Interleukin-33 (IL-33), a member of the IL-1 family of cytokines, is expressed by many cell types following pro-inflammatory stimulation and is thought to be released on cell lysis. The 30 kDa human IL-33 is converted by CASP1 to a 18 kDa protein. IL-33 binds to and signals through ST2 (IL1R1) and its stimulation recruits MYD88, IRAK, IRAK4, and TRAF6, followed by phosphorylation of ERK1 (MAPK3)/ERK2 (MAPK1), p38 (MAPK14), and JNK. The ability of IL-33 to target numerous immune cell types, like Th2-like cells, mast cells, and B1 cells, and to induce cytokine and chemokine production underlines its potential in influencing the outcome of a wide range of diseases, such as arthritis, asthma, atopic allergy and anaphylaxis, cardiovascular disease/atherosclerosis, nervous system diseases, and sepsis.



Lane 1: IL-33 (human recombinant; His-tagged)
Lane 2: IL-33 (human recombinant; FLAG®-tagged)

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10809

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/17/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com