

# PRODUCT INFORMATION



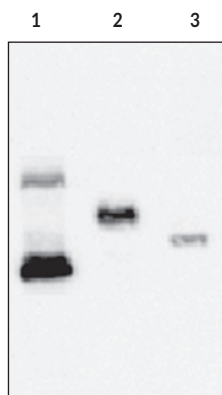
## Interleukin-33 Monoclonal Antibody (Clone IL333068A)

Item No. 10810

### Overview and Properties

|                          |                                                                                                                                                                                                                                                 |
|--------------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Contents:</b>         | This vial contains 50 or 100 µg of protein G-purified antibody.                                                                                                                                                                                 |
| <b>Synonyms:</b>         | IL-1F11, IL-33, NF-HEV                                                                                                                                                                                                                          |
| <b>Immunogen:</b>        | Recombinant human IL-33                                                                                                                                                                                                                         |
| <b>Cross Reactivity:</b> | (+) Human IL-33, murine IL-33 (weakly)                                                                                                                                                                                                          |
| <b>Form:</b>             | Liquid                                                                                                                                                                                                                                          |
| <b>Storage:</b>          | -20°C (as supplied)                                                                                                                                                                                                                             |
| <b>Stability:</b>        | ≥1 year                                                                                                                                                                                                                                         |
| <b>Storage Buffer:</b>   | PBS, pH 7.4                                                                                                                                                                                                                                     |
| <b>Clone:</b>            | IL333068A                                                                                                                                                                                                                                       |
| <b>Host:</b>             | Mouse                                                                                                                                                                                                                                           |
| <b>Isotype:</b>          | IgG2bk                                                                                                                                                                                                                                          |
| <b>Applications:</b>     | ELISA and Western blot (WB); the recommended starting dilution for ELISA (direct or indirect) and WB is 1:2,000-1:5,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

### Image



Lane 1: Recombinant IL-33 (human) (His-tagged)  
Lane 2: Recombinant IL-33 (human) (FLAG<sup>®</sup>-tagged)  
Lane 3: Recombinant IL-33 (mouse) (FLAG<sup>®</sup>-tagged)

### Description

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 B<sub>1</sub> 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.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/10/2022

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM