

PRODUCT INFORMATION



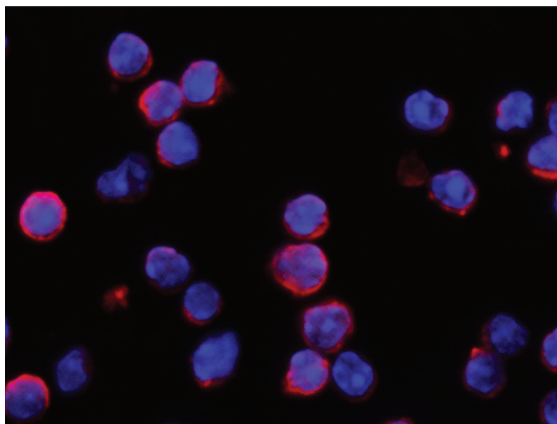
IFNAR1 Rabbit Monoclonal Antibody

Item No. 38087

Overview and Properties

| | |
|----------------------------|--|
| Contents: | This vial contains 50 or 100 µl of protein A-affinity purified monoclonal antibody. |
| Synonyms: | Interferon Alpha/Beta Receptor 1, Interferon α/β Receptor 1, CRF2-1, CRF2 member 1, Cytokine Receptor Class-II Member 1, Cytokine Receptor Family 2 Member 1, Type I IFN Receptor 1, Type I Interferon Receptor 1 |
| Immunogen: | Recombinant mouse IFNAR1 |
| Cross Reactivity: | (+) <i>Ifnar1</i> |
| Species Reactivity: | (+) Mouse; other species not tested |
| Form: | Liquid |
| Storage: | 80°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | 0.2 µm filtered solution in PBS |
| Clone: | 110 |
| Host: | Rabbit |
| Isotype: | IgG |
| Applications: | Immunocytochemistry (ICC) and Immunofluorescence (IF); the recommended starting concentration is 1:20-1:100 for ICC and IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Image



Immunofluorescent analysis of mouse IFNAR1 in mouse splenocytes. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with IFNAR1 Rabbit Monoclonal Antibody at a dilution of 1:60 at 37°C for one hour. Cells were then stained with an Alexa Fluor® 594-conjugated goat anti-rabbit IgG secondary antibody (red) and counterstained with DAPI (blue).

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, 02/08/2023

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

PRODUCT INFORMATION



Description

α/β receptor 1 (IFNAR1), also known as type I IFN receptor 1, is a class II helical cytokine receptor involved in the antiviral response.¹ It is composed of an extracellular domain, which contains four fibronectin type III subdomains, a transmembrane domain, and an intracellular domain, which contains a tyrosine kinase 2 (TYK2) phosphorylation site, and forms a heterodimer with IFNAR2.² It is ubiquitously expressed and binds to type I IFNs, including IFN- α , IFN- β , and IFN- ω , which, upon binding, together with IFNAR2 signals through JAK1, TYK2, and various STATs.³ Knockout of *Ifnar1* sensitizes fibroblasts to severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) infection *in vitro* and has commonly been used as an immunosuppressive model to study viral infections *in vivo*.^{4,5} SNPs in *IFNAR1* are associated with increased risk of viral infections, lung cancer, multiple sclerosis, and cerebral malaria, as well as adverse reactions to vaccinations.¹ Cayman's IFNAR1 Rabbit Monoclonal Antibody can be used for immunocytochemistry (ICC) and immunofluorescence (IF) applications.

References

1. de Weerd, N.A., J.P., V., Lim, S.S., *et al.* Structural integrity with functional plasticity: What type I IFN receptor polymorphisms reveal. *J. Leukoc. Biol.* **108(3)**, 909-924 (2020).
2. Borden, E.C., Sen, G.C., Uze, G., *et al.* Interferons at age 50: Past, current and future impact on biomedicine. *Nat. Rev. Drug Discov.* **6(12)**, 975-990 (2007).
3. de Weerd, N.A. and Nguyen, T. The interferons and their receptors - distribution and regulation. *Immunol. Cell Biol.* **90(5)**, 483-491 (2012).
4. Zhang, Q., Bastard, P., Liu, Z., *et al.* Inborn errors of type I IFN immunity in patients with life-threatening COVID-19. *Science* **370(6515)**, eabd4570 (2020).
5. Marín-Lopez, A., Calvo-Pinilla, E., Moreno, S., *et al.* Modeling arboviral infection in mice lacking the interferon alpha/beta receptor. *Viruses* **11(1)**, 35 (2019).