

IFNGR1 recombinant monoclonal antibody (062)

This antibody is used in research focusing on immunological pathways, cytokine signaling, and immune regulation.

Interferon- γ receptor 1 (IFNGR1) is a key component of the receptor complex for interferon-gamma (IFN- γ), which is critical for immune responses against viral and bacterial infections. This antibody is used in research focusing on immunological pathways, cytokine signaling, and immune regulation.

This antibody is covered by our [Worry-Free Guarantee](#).

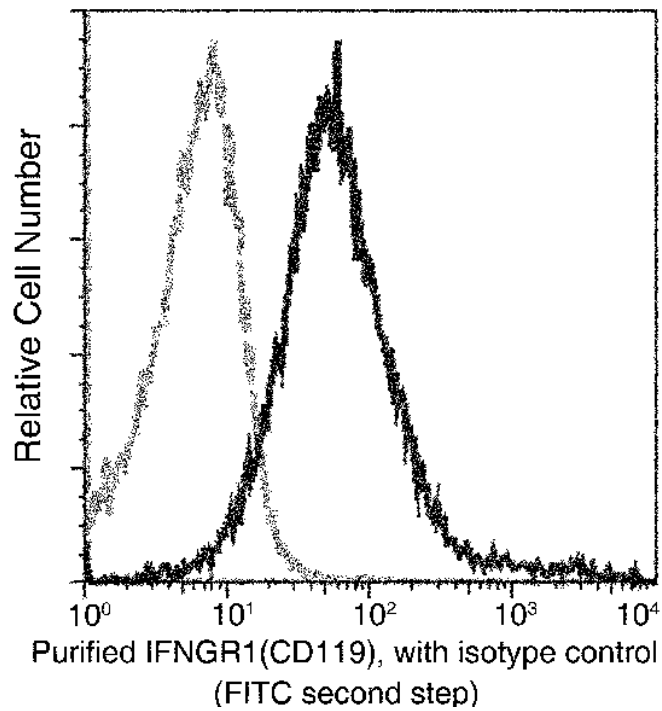
Ordering Information

[Order Online »](#)

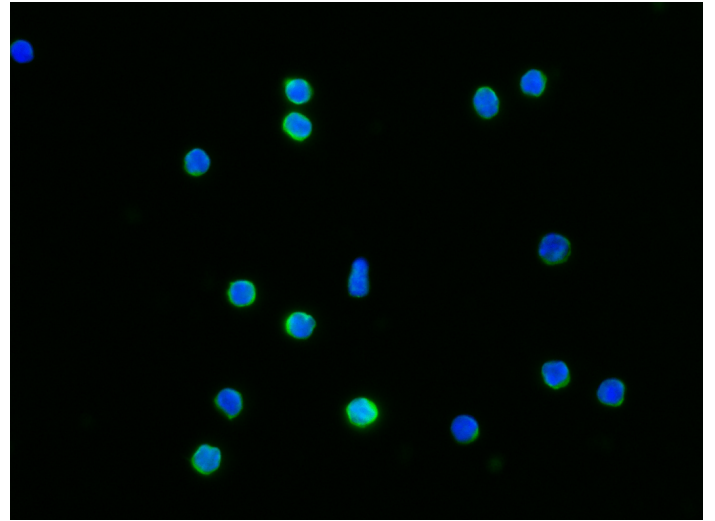
ENZ-ABS1033-0100	100 μ l
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse IFNGR1(CD119) expression on BABL/c splenocytes. Cells were stained with purified anti-Mouse IFNGR1(CD119), then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of mouse IFNGR1 in mouse splenocytes. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-mouse IFNGR1 monoclonal antibody (1:60) at 4°C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue).

Handling & Storage

Handling	Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	IFN-γ receptor, Interferon-γ receptor 1, CD119
Application	ELISA, Flow Cytometry, ICC, IF
Clone	062
Crossreactivity	No cross-reactivity in ELISA with human LDLR.
Formulation	Liquid. 0.2 μm filtered solution in PBS. Preservative free. Carrier-free.
Host	Rabbit
Immunogen	Recombinant mouse IFNGR1.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunocytochemistry / Immunofluorescence (1:20-1:100)Flow Cytometry (1:25-1:100)ELISA (Capture) (1:250-1:2'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: October 7, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)