

IL-1R3 monoclonal antibody (7E4G1E8)

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

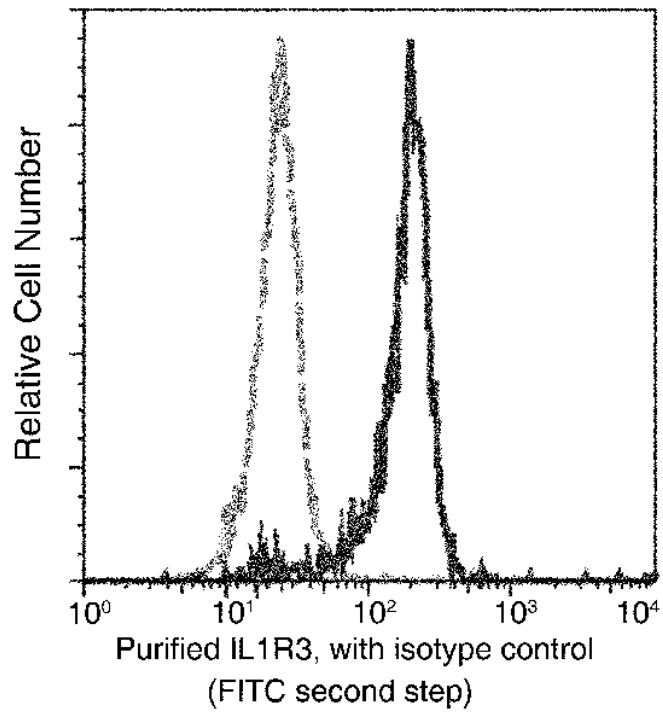
Ordering Information

[Order Online »](#)

ENZ-ABS935-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human IL1R3 expression on human whole blood monocytes. Cells were stained with purified anti-Human IL1R3, then a FITC-conjugated second step antibody. The histogram were derived from gated events with the forward and side light-scatter characteristics of viable monocytes.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Interleukin-1 receptor 3, IL-1RAcP, Interleukin-1 receptor accessory protein
Application	ELISA, Flow Cytometry
Clone	7E4G1E8
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Mouse
Immunogen	Recombinant human IL1R3 protein.
Isotype	IgG2a
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	ELISA (1:1'000-1:2'000)Flow Cytometry(1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: September 17, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430
info-
usa@enzolifesciences.com

European Sales Office

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands

& Luxembourg
Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France

Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany

Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland

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