CD213A2 recombinant monoclonal antibody (018)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

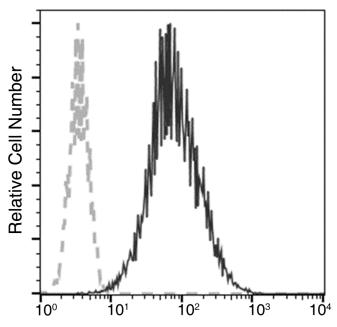
Order Online »

ENZ-ABS943-0100

100µl

Manuals, SDS & CofA

View Online »



Purified IL13Ra2(CD213a2), with isotype control (FITC second step)

Flow cytometric analysis of Human IL13Ra2(CD213a2) expression on A375 cells. Cells were stained with purified anti-Human IL13Ra2(CD213a2), 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.

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 IL13RA2, Interleukin-13 receptor subunit α -2

Application Flow Cytometry

018 Clone

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human IL13RA2 / CD213A2 protein.

IgG Isotype

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions 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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

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

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