

Mouse IgG1 isotype control, monoclonal antibody (MOPC-21) (FITC conjugate)

Isotype matched controls are useful in flow cytometry to determine a gross estimate of non-specific binding that may be observed with a specific antibody of interest. In general, an isotype control should be of the same species and isotype as the antibody of interest, used at the same concentration, and conjugated to the same fluorophore. This reagent is an IgG1 isotype mouse monoclonal antibody (Clone # MOPC-21), conjugated to FITC. It is intended for use as an isotype control antibody for flow cytometry.

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

Citations: 20

[View Online »](#)

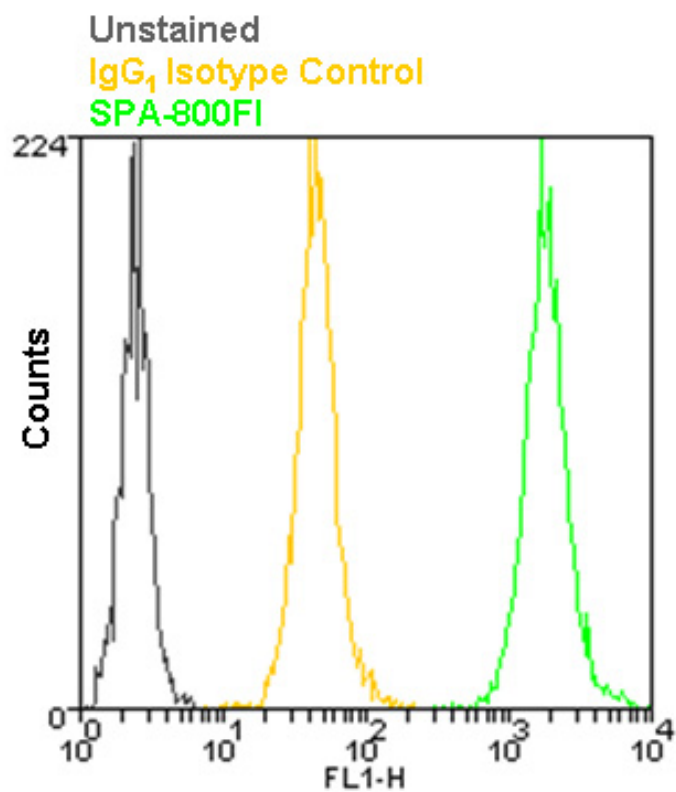
Ordering Information

[Order Online »](#)

ADI-SAB-600FI-050	50µg
ADI-SAB-600FI-200	200µg

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10^6 Jurkat cells using the Isotype Control for IgG1 FITC (Prod. No. ADI-SAB-600FI) compared with Hsp27, mAb (Prod. No. ADI-SPA-800FI), at a concentration of $50\mu\text{g/ml}$.

Handling & Storage

Handling Avoid freeze/thaw cycles. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry

Clone MOPC-21

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Host Mouse

Isotype IgG1

Purity Detail Protein G affinity purified.

Source Purified from ascites.

Species Reactivity Species independent

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
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