

# Mouse IgG<sub>1</sub> isotype control, monoclonal antibody (MOPC-21)

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). It is intended for use as an isotype control antibody for flow cytometry.

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

Citations: 1 [View Online »](#)

Ordering Information [Order Online »](#)

ADI-SAB-600-0050	50µg
ADI-SAB-600-0200	200µg

Manuals, SDS & CofA [View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°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.

**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](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)