

# Mouse IgG<sub>1</sub> isotype control, monoclonal antibody (MOPC-21) (DyLight™ 488 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 DyLight™-488. It is intended for use as an isotype control antibody for flow cytometry.

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

Citations: 2

[View Online »](#)

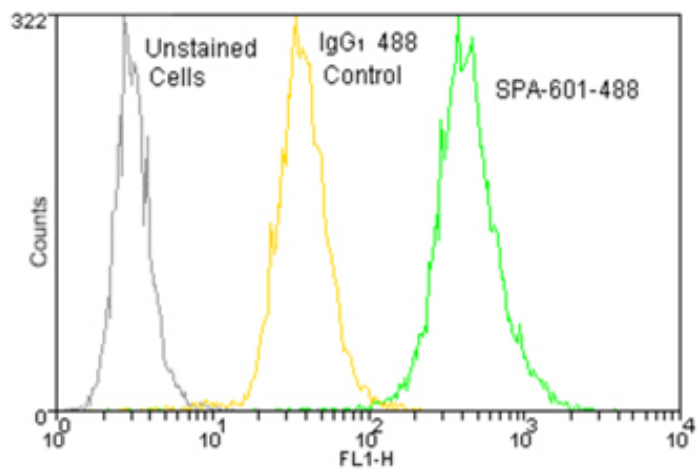
## Ordering Information

[Order Online »](#)

ADI-SAB-600-488-D	50µg
ADI-SAB-600-488-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10<sup>6</sup> Jurkat cells using the Isotype Control for IgG1 488 (Prod. No. ADI-SAB-600-488) compared with a Calreticulin, mAb (Prod. No. ADI-SPA-601-488), at a concentration of 50µg/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](#).



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-  
fr@enzolifesciences.com](mailto:info-fr@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)