

Human IgG1 Isotype Control

Human IgG1 are purified from human plasma by fractionation, ion exchange and affinity chromatography. The purified IgG represent a single subclass and light chain type (κ). Source material has been shown to be non reactive for HBsAg, anti-HCV, anti-HBc, and negative for anti-HIV 1 & 2 by FDA approved tests. Negative control in experiments with recombinant proteins fused to human IgG1 Fc.

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

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-133-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored as recommended.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, Flow Cytometry, IP, WB
Application Notes	Negative control in assays where receptors:Fc (IgG1) are used.
Formulation	Lyophilized from PBS, pH 7.2.
Host	Human
Isotype	IgG1
Purity Detail	Purified from myeloma plasma.
Reconstitution	Reconstitute with 100µl sterile water.
Source	Purified from human plasma.
Species Reactivity	Species independent
Technical Info / Product Notes	Cited samples: For an overview on cited samples please click here.
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-
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)