

Mouse Control IgG1

Isotype control for mouse IgG1 antibodies
Catalog code: mabg1-ctrlm, mabg1-ctrlm-05
<https://www.invivogen.com/control-igg1>

For research use only
Version 23L12-MM

PRODUCT INFORMATION

Contents: Mouse Control IgG1 purified monoclonal antibody (mAb) is provided azide-free and lyophilized. It is available in two quantities:

- mabg1-ctrlm:** 100 µg Mouse Control IgG1
- mabg1-ctrlm-05:** 5 x 100 µg Mouse Control IgG1

Target: *E. coli* β-galactosidase (β-Gal)

Species reactivity: Does not react with murine proteins

Clone: T8E5

Isotype: Mouse IgG1

Source: Hybridoma

Light chain type: Kappa

Purification: By affinity chromatography

Formulation: 0.2 µm filtered solution in a sodium phosphate buffer with glycerine, saccharose, and stabilizing agents

Tested applications: Mouse Control IgG1 is provided as a negative control for use in conjunction with InvivoGen's mouse IgG1 monoclonal antibodies.

Antibody resuspension (0.1 mg/ml)

Note: Ensure you see the lyophilized pellet before resuspension.

Resuspend Mouse Control IgG1 with sterile water:

Add 1 ml of sterile water per 100 µg vial.

Storage and stability

- Product is shipped at room temperature. Store lyophilized antibody at -20 °C.
- Reconstituted antibody is stable for 1 month when stored at 4°C and for 1 year when aliquoted and stored at -20 °C. Avoid repeated freeze-thaw cycles.

Quality control

- Functionality of Mouse Control IgG1 has been tested using flow cytometry.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Mouse Control IgG1 is a monoclonal mouse antibody specific for *E. coli* β-galactosidase (β-Gal). This antibody was obtained from a mouse immunized with DNA encoding β-Gal. This monoclonal antibody was purified by affinity chromatography.

APPLICATIONS

Mouse Control IgG1 is provided as a negative control for use in conjunction with InvivoGen's mouse IgG1 monoclonal antibodies. This isotype control product should be used at the same concentration as the specific monoclonal antibody. Optimal dilutions should be determined for each application.

RELATED PRODUCTS

Product	Cat. Code
Mouse Control IgG2a	mabg2a-ctrlm

InvivoGen provides a collection of monoclonal antibodies targeting human and murine TLRs (Toll-Like receptors) and cytokines. For more information, visit <https://www.invivogen.com/antibodies-list>.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com

