

Mouse IgG2b Control

Isotype control for mouse IgG2b antibodies

Catalog # mabg2b-ctrlm

For research use only, not for diagnostic or therapeutic use

Version # 15C23-MM

PRODUCT INFORMATION

Content

100 μ g purified Mouse IgG2b Control antibody, provided azide-free and lyophilized

Specificity: β -galactosidase

Clone: T8H6

Isotype: Mouse IgG2b

Formulation: 0.2 μ m filtered solution in phosphate/glycine with 5% saccharose

Antibody resuspension

Add 1 ml of sterile water to obtain a concentration of 0.1 mg/ml.

Storage

- Product is shipped at room temperature. Store lyophilized antibody at -20 °C. Product is stable for 6 months.

- Reconstituted antibody is stable 1 month when stored at 4 °C and 6 months when aliquoted and stored at -20 °C. Avoid repeated freeze-thaw cycles.

Quality control

- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) is confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Mouse IgG2b Control 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.

APPLICATION

Mouse IgG2b Control is provided as a negative control for use in conjunction with InvivoGen's mouse IgG2b 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. This antibody should not be used at a concentration higher than 500 ng/10⁶ cells when used as a negative control for flow cytometry.

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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com

