

Rat IgG2b Control

Isotype Control for Rat IgG2b

Catalog # mabg2b-ctrlt

For research use only, not for diagnostic or therapeutic use

Version # 14B27-MM

PRODUCT INFORMATION

Content

100 µg purified Rat IgG2b Control antibody, provided azide-free and lyophilized

Clone: C1-12G10

Isotype: Rat IgG2b

Formulation: 0.2 µm 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.

Description

Rat IgG2b Control is a monoclonal rat antibody specific for *E. coli* β-galactosidase (β-Gal). This antibody was obtained from a rat immunized with DNA encoding β-Gal. This monoclonal antibody was purified by affinity chromatography.

Specificity

This antibody was purified for use as a rat IgG2b isotype control. This antibodies recognizes β-Gal.

APPLICATION

Rat IgG2b Control is provided as a negative control for use in conjunction with InvivoGen's rat 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.

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



www.invivogen.com