

Rat IgG2a Control

Isotype Control for Rat IgG2a

Catalog # mabg2a-ctrlt

For research use only, not for diagnostic or therapeutic use

Version # 14B27-MM

PRODUCT INFORMATION

Content

100 μ g purified Rat IgG2a Control antibody, provided azide-free and lyophilized

Clone: C1-2A7

Isotype: Rat IgG2a

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.

Description

Rat IgG2a 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 IgG2a isotype control. This antibodies recognizes β -Gal.

APPLICATION

Rat IgG2a Control is provided as a negative control for use in conjunction with InvivoGen's rat IgG2a 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

