

Control-IgA2

Human IgA2 Isotype Control

Catalog # maba2-ctrl

For research use only, not for diagnostic or therapeutic use

Version # 09G01-MT

PRODUCT INFORMATION

Content

100 μ g purified Control-IgA2 antibody, provided azide-free and lyophilized

Clone: T9C6

Isotype: Human IgA2

Formulation: 0.2 μ m filtered solution in Tris/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

Control-IgA2 is a chimeric monoclonal antibody specific for *E. coli* β -galactosidase (β -Gal). It was generated by combining the constant domains of the human IgA2 molecule with the murine variable regions of a monoclonal IgG antibody obtained from a mouse immunized with DNA encoding β -Gal. This monoclonal antibody was purified by affinity chromatography.

Specificity

This antibody was purified for use as a human IgA2 isotype control. This antibodies recognizes β -Gal.

APPLICATION

Control-IgA2 is provided as a negative control for use in conjunction with InvivoGen's IgA2 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

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite A
San Diego, CA 92121 - USA