

# Human IgA2 Control

Isotype control for human IgA2 antibodies

Catalog # maba2-ctrl

For research use only, not for diagnostic or therapeutic use

Version # 15C23-MM

## PRODUCT INFORMATION

### Content

100  $\mu$ g purified Human IgA2 control antibody, provided azide-free and lyophilized

**Specificity:**  $\beta$ -galactosidase

**Clone:** T9C6

**Isotype:** Human IgA2

**Formulation:** 0.2  $\mu$ 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  $-20^{\circ}\text{C}$ . Product is stable for 6 months.

- Reconstituted antibody is stable 1 month when stored at  $4^{\circ}\text{C}$  and 6 months when aliquoted and stored at  $-20^{\circ}\text{C}$ . 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

Human IgA2 Control is a chimeric monoclonal antibody specific for *E. coli*  $\beta$ -galactosidase ( $\beta$ -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  $\beta$ -Gal. This monoclonal antibody was purified by affinity chromatography.

## APPLICATION

Human IgA2 Control 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](mailto:info@invivogen.com)

Website: [www.invivogen.com](http://www.invivogen.com)



3950 Sorrento Valley Blvd. Suite 100  
San Diego, CA 92121 - USA