

# Human IgA2 Control

Isotype control for human IgA2 antibodies

Catalog code: maba2-ctrl, maba2-ctrl-05

<https://www.invivogen.com/control-iga>

For research use only

Version 23L15-MM

## PRODUCT INFORMATION

**Contents:** Human IgA2 Control purified monoclonal antibody (mAb) is provided azide-free and lyophilized. It is available in two quantities:

**maba2-ctrl:** 100 µg Human IgA2 Control

**maba2-ctrl-05:** 5 x 100 µg Human IgA2 Control

**Target:** *E. coli* β-galactosidase (β-Gal)

**Species reactivity:** Does not react with human proteins

**Clone:** T9C6

**Isotype:** Human IgA2

**Source:** CHO cells

**Light chain type:** Kappa

**Purification:** By affinity chromatography

**Formulation:** 0.2 µm filtered solution in TRIS with glycine, saccharose, and stabilizing agents

**Tested applications:** Human IgA2 Control is provided as a negative control for use in conjunction with InvivoGen's human IgA2 monoclonal antibodies.

### Antibody resuspension (0.1 mg/ml)

*Note:* Ensure you see the lyophilized pellet before resuspension.

Resuspend Human IgA2 Control with sterile water:

Add 1 ml of sterile water per 100 µg vial.

### Storage and stability

- Product is shipped at room temperature. Store lyophilized antibody at -20°C.

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

### Quality control

- Functionality of Human IgA2 Control has been tested using flow cytometry.

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

## DESCRIPTION

Human IgA2 Control is a mono-glycosylated monoclonal human antibody specific for *E. coli* β-galactosidase (β-Gal). This antibody features the constant region of the human IgA2 isotype and the variable region of Mouse IgG2a Control (clone: T9C6) targeting *E. coli* β-Gal. Therefore, it does not react with human targets. This monoclonal antibody was purified by affinity chromatography.

## APPLICATIONS

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

## RELATED PRODUCTS

Product	Cat. Code
Anti-hTLR2-IgA	maba2-htlr2-2
Anti-hTLR5-IgA	maba2-htlr5-2

InvivoGen provides a collection of monoclonal antibodies targeting human and murine TLRs (Toll-Like receptors) and cytokines. For more information, visit <https://www.invivogen.com/antibodies-list>.

## 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 Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

