

Anti-β-Gal-hIgG4

Human IgG4 monoclonal antibody against β-galactosidase - Isotype control

Catalog # bgal-mab4

For research use only, not for diagnostic or therapeutic use

Version # 16E03-MM

PRODUCT INFORMATION

Content: 200 µg Anti-β-Gal-hIgG4, purified monoclonal antibody, provided azide-free and lyophilized

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

Isotype: Human IgG4

Source: CHO cells

Formulation: 0.2 µm filtered solution in sodium phosphate buffer with glycine, saccharose and stabilizing agents.

Antibody resuspension

Add 2 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. Lyophilized product is stable for 1 year.

- 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

- Absence of binding of Anti-β-Gal-hIgG4 to human cell lines has been tested using flow cytometry.

- The complete sequence of this antibody has been verified.

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

DESCRIPTION

Anti-β-Gal-hIgG4 features the constant region of the human IgG4 isotype and the variable region of [Mouse IgG2a Control](#). Mouse IgG2a Control is a mouse IgG2a monoclonal antibody that targets *E. coli* β-galactosidase (β-Gal). This antibody was generated by DNA immunization with a plasmid expressing the β-Gal gene in Swiss mice. It can be used as an isotype control for human IgG4 antibodies.

Human IgG4 is the least common IgG present in serum. Human IgG4 displays low antibody-dependent cell-mediated cytotoxicity (ADCC) and no complement-dependent cytotoxicity (CDC).

Anti-β-Gal-hIgG4 was generated by recombinant DNA technology. It has been produced in CHO cells and purified by affinity chromatography.

ANTIBODY ISOTYPE FAMILY

For your research, InvivoGen provides an anti-β-Gal isotype family. This family consists of monoclonal antibodies comprising the variable region of a mouse monoclonal antibody targeting *E. coli* β-galactosidase (β-Gal), named mouse IgG2a control, and the constant region of three different human isotypes; IgG1, IgG4 and IgA2. The isotypes differ in their functional locations and effector functions, such as complement-dependent cytotoxicity (CDC) and antibody-dependent cell-mediated cytotoxicity (ADCC), as presented in the table below.

Isotype	Description
Human IgG1	Most abundant IgG present in serum High CDC, high ADCC
Human IgG4	Least common IgG present in serum No CDC, low ADCC
Human IgA2	Major class in secretions, oligomeric forms, highly resistant to enzymatic degradation. No CDC, low ADCC

RELATED PRODUCTS

Product	Catalog Code
Anti-β-Gal-hIgG1	bgal-mab1
Anti-β-Gal-hIgA2	hbgal-mab7
Mouse IgG2a Control	mabg2a-ctrlm

Several antibody isotype families are available, such as Anti-hCD20, Anti-HER2 and Anti-hPD1.

For more information visit www.invivogen.com/antibody-isotypes

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 3-622-34-80

E-mail: info@invivogen.com