

Anti- β -Gal-hIgG1*

Recombinant human monoclonal IgG1 antibody against β -galactosidase

Catalog code: bgal-mab1-1

<https://www.invivogen.com/anti-bgal-0>

For research use only

Version 23J12-NJ

PRODUCT INFORMATION

Contents: 1 mg purified Anti- β -Gal-hIgG1 monoclonal antibody (mAb), provided azide-free and lyophilized

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

Source: Chinese hamster ovary (CHO) cells

Isotype: Human IgG1, kappa

Purification: By affinity chromatography with protein A

Formulation: 0.2 μ m filtered solution in 150 mM sodium chloride, 20 mM sodium phosphate buffer with 5% saccharose

Tested applications: Flow cytometry, ELISA, Antibody-drug conjugation

Storage

- Product is shipped at room temperature. Store lyophilized antibody at -20 °C. Lyophilized product is stable for at least 1 year.
- Reconstituted antibody is stable for 1 month at 4 °C and for 1 year at -20 °C. Avoid repeated freeze-thaw cycles.

Quality control

- Binding of Anti- β -Gal-hIgG1* to β -galactosidase has been validated using ELISA.
- The absence of non-specific binding of Anti- β -Gal-hIgG1* has been validated using flow cytometry and antibody-drug conjugate (ADC)-based cellular assays.
- The complete sequence of the antibody has been verified.
- Absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and TLR4 cellular assays.

PRODUCT DESCRIPTION

Anti- β -Gal-hIgG1* is a recombinant monoclonal antibody (mAb) featuring the variable region of a murine mAb which targets *E. coli* β -galactosidase, and the constant region of the hIgG1 isotype. Anti- β -Gal-hIgG1* was generated by recombinant DNA technology, produced in CHO cells, and purified by affinity chromatography with protein A.

IgG1 Isotype effector function

Human IgG1 binds with high affinity to the Fc receptor on phagocytic cells. Therefore, hIgG1 mAbs display high effector function, including antibody-dependent cell-mediated cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC).

APPLICATIONS

Anti- β -Gal-hIgG1* has been specifically developed for the generation of antibody-drug conjugates (ADCs). The antibody formulation contains no tris, cysteine, glycine, nor any other amine or thiol function.

This antibody can be used to generate control ADCs which do not recognize any antigen other than β -galactosidase.

For examples of ADC-mediated cell based assays, please refer to the technical and validation documents of InvivoGen's STING and TLR7 agonists for bioconjugation, as well as Anti-HER2-hIgG1 and Anti-TROP2-hIgG1 mAbs for ADC:

<https://www.invivogen.com/sting-conjugatable-ligands>

<https://www.invivogen.com/tlr7-conjugatable-ligands>

<https://www.invivogen.com/anti-her2-0>

<https://www.invivogen.com/anti-trop2>

To learn more about antibody drug conjugates and conjugatable ligands, please visit <https://www.invivogen.com/conjugatable-ligands>.

RELATED PRODUCTS

Product	Description	Cat.Code
TL7-887	Conjugatable TLR7 ligand	vac-tl7887
TL7-975	Conjugatable TLR7 ligand	vac-tl7975
STG-982	Conjugatable STING ligand	vac-stg982
STG-968	Conjugatable STING ligand	vac-stg968
Anti-HER2-hIgG1	Monoclonal antibody	her2-mab1-1
Anti-TROP2-hIgG1	Monoclonal antibody	trop2-mab1-1

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