

# Anti-TROP2-hIgG1

Recombinant human monoclonal IgG1 antibody against TROP2

Catalog code: trop2-mab1-1

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

For research use only

Version 23J12-NJ

## PRODUCT INFORMATION

**Contents:** 1 mg purified Anti-TROP2-hIgG1 monoclonal antibody (mAb), provided azide-free and lyophilized

**Target:** Human trophoblast cell-surface antigen 2 (TROP2)

**Variable region biosimilar:** Sacituzumab

**Source:** Chinese hamster ovary (CHO) cells

**Isotype:** Human IgG1, kappa

**Purification:** By affinity chromatography with protein A

**Formulation:** 0.2 µm 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-TROP2-hIgG1 to TROP2 has been validated using flow cytometry, ELISA, 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-TROP2-hIgG1 is a recombinant monoclonal antibody (mAb) that targets human trophoblast cell-surface antigen 2 (TROP2, also known as tumor associated calcium signal transducer 2, TACSTD2). It features the variable region of sacituzumab and the constant region of the hIgG1 isotype. Anti-TROP2-hIgG1 was generated by recombinant DNA technology, produced in CHO cells, and purified by affinity chromatography with protein A.

### IgG1 Isotype effector function

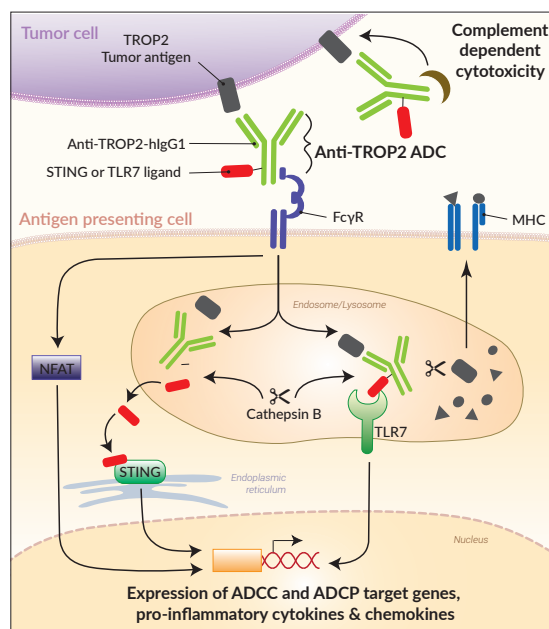
Human IgG1 binds with high affinity to the Fc receptor on phagocytic cells. Therefore, hIgG1 mAbs display high effector function, including antibody-dependent cell-mediated cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC).

### Applications

Anti-TROP2-hIgG1 has been specifically developed for the generation of antibody-drug conjugates. The antibody formulation contains no tris, cysteine, glycine, nor any other amine or thiol function.

## BACKGROUND

TROP2 (human trophoblast cell-surface antigen 2) is a transmembrane protein expressed in normal human tissues, including breast, cervix, kidney, lung, and pancreas, and is upregulated in most tumors of epithelial origin<sup>1</sup>. TROP2 plays an important role in embryonic development and is implicated in several oncogenic signaling pathways leading to proliferation<sup>1</sup>. One therapeutic strategy aiming at killing TROP2<sup>+</sup> cancer cells, notably in triple-negative breast cancer patients, uses the FDA-approved antibody-drug conjugate (ADC) Trodelvy<sup>2,3</sup>. It consists of the combination of the Anti-TROP2 sacituzumab and the SN-38 cytotoxic payload<sup>2,3</sup>. This ADC allows targeted drug delivery to cancer cells (thus limiting systemic side-effects) in addition to the mAb-mediated effector functions<sup>2,3</sup>. Other TROP2-targeting ADCs featuring different cytotoxic cargos have been developed and have progressed to human clinical trials<sup>4</sup>.



1. Trerotola, M., et al., 2013. Upregulation of Trop-2 quantitatively stimulates human cancer growth. *Oncogene* 32, 222.
2. Drago J.Z. et al., 2021. Unlocking the potential of antibody-drug conjugates for cancer therapy. *Nat Rev Clin Oncol*. 18-6:327.
3. Tarantino P. et al., 2022. Antibody-drug conjugates: smart chemotherapy delivery across tumor histologies. *CA Cancer J Clin*. 72:165.
4. Shaffer, C. 2021. Trop2 deal heats up antibody-drug conjugate space in cancer. *Nat Biotechnol* 39, 128.

## 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)

## PROTOCOLS

### Antibody-drug conjugation

For conjugation protocols of the monoclonal antibody to STING or TLR7 agonists, using either cysteine-based thioether coupling or click-chemistry coupling, please refer to the technical data sheets of :

- STG-982 and STG-968, STING agonists with maleimide and azido group, respectively: <https://www.invivogen.com/sting-conjugatable-ligands>.
- TL7-887 and TL7-975, TLR7 agonists with maleimide and azido group, respectively: <https://www.invivogen.com/tlr7-conjugatable-ligands>.

*Note: Anti- $\beta$ -Gal-hIgG1\* mAb can be used to generate control ADC.*

### ADC-mediated cell stimulation

The biological functions of ADCs can be assessed using reporter cell lines co-expressing the targeted tumor antigen and PRR (e.g. STING or TLR7). Examples of such mono-culture models are shown in the validation data sheets of the STG-968 and TL7-975 conjugatable ligands (see URLs above).

Alternatively, the biological functions of ADCs can be assessed using co-cultures of tumor cell lines expressing the targeted tumor antigen and human peripheral blood monocytes (PBMCs), myeloid antigen presenting cells, or reporter THP-1 monocytes. Examples of such co-culture models are shown in the validation data sheets of the Anti-TROP2-hIgG1, as well as in different reports<sup>6,7</sup>.

6. Ackerman S.E. *et al.*, 2021. Immune-stimulating antibody conjugates elicit robust myeloid activation and durable antitumor immunity. *Nature Cancer*. 2(1):18-33.

7. Duvall J.R. *et al.*, 2021. XMT-2056, a well-tolerated, Immunosynthen-based STING-agonist antibody-drug conjugate which induces anti-tumor immune activity. *Cancer Research* 81(13):1738.

## 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- $\beta$ -Gal-hIgG1*	Monoclonal antibody	bgal-mab1-1
Anti-HER2-hIgG1	Monoclonal antibody	her2-mab1-1
HEK-Blue™ IL-6 cells	IL-6 reporter cells	hkb-hil6
QUANTI-Blue™	Detection reagent	rep-qbs

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

### 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)