

bsAb CD3-CD28 for T cell expansion

Bispecific antibody (bsAb) against human CD3 and human CD28

Catalog codes: bsab-tex-1, bsab-tex-2, bsab-tex-4

For research use only

Version 24J01-NJ

PRODUCT INFORMATION

Contents: bsAb CD3-CD28 is a purified bispecific antibody (bsAb), provided azide-free and lyophilized. It is available in three quantities:

- **bsab-tex-1:** 100 µg
- **bsab-tex-2:** 2 x 100 µg
- **bsab-tex-4:** 4 x 100 µg

Target: Human (h)CD3 (clone OKT3) and hCD28 (clone 15E8)

Format: IgG-like (scFv)₂-(scFv)₂-Fc molecule

Source: Chinese hamster ovary (CHO) cells

Capacity: 100 µg of bsAb CD3-CD28 is sufficient to activate and expand up to 1x10⁸ enriched T cells or peripheral blood mononuclear cells (PBMCs), when used at the recommended titer of 1:100.

Storage and stability

- Product is shipped at room temperature. Upon receipt, store lyophilized antibody at -20 °C. Lyophilized product is stable for at least 1 year.
- Reconstituted antibody is stable for 1 year when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.

Quality control

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

Antibody resuspension (0.1 mg/ml)

Note: Ensure you see the lyophilized pellet before resuspension.

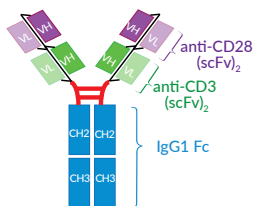
- Add 1 ml of sterile H₂O to the vial and gently pipette until completely resuspended.

APPLICATION

The bsAb CD3-CD28 is a bispecific antibody (bsAb) designed to activate and expand enriched T cell populations or resting T cells from PBMCs. T cell expansion is achieved by culturing for up to 1-2 months.

PRODUCT INFORMATION

The bispecific antibody (bsAb) CD3-CD28 is a fusion protein dimer comprising tandem single-chain variable fragments (scFv)₂ from two monoclonal antibodies (mAbs). These mAbs were derived from the clones OKT3 and 15E8 targeting the human CD3 and CD28 receptors, respectively. This so-called (scFv)₂-(scFv)₂-Fc molecule belongs to the family of IgG-like antibodies. It features an IgG1 Fc fragment, which contributes to increased solubility, serum half-life, as well as facilitated purification.



CELLULAR ASSAY

All steps in the protocol have to be performed under sterile conditions. *Note: For optimal T cell expansion, a daily inspection of culture is required.*

Required Cell Culture Medium

- **Growth Medium:** RPMI, 10% (v/v) heat-inactivated FBS, 100 U/ml penicillin, 100 µg/ml streptomycin, 50 µM β-mercaptoethanol, 1 mM Sodium pyruvate, 1x NEAA

Required Conditions

- 5% CO₂, 37°C Incubator

Required Supplements

- [Recombinant human IL-2 cytokine](#)

Day 1: T cell isolation and activation

1. Isolate T cells according to your protocol.
2. Determine cell number.
3. Resuspend T cells at 1x10⁶ cells/ml in freshly prewarmed media (*see table on page 2*).
4. Add [recombinant human IL-2](#) (final concentration 20 IU/ml) and bsAb CD3-CD28 (final concentration **1 µg/ml**) directly to the cells.
5. Incubate at 37 °C, 5% CO₂ for **2 days**.

Day 3: Stop activation

1. Remove residual reagent bsAb CD3-CD28 by centrifugation at 300 x g (RCF) for 5 min.
2. Aspirate supernatant.
3. Determine cell number.
4. Resuspend T cells at 1x10⁶ cells/ml in freshly prewarmed medium.
5. Add [recombinant human IL-2](#) (final concentration 20 IU/ml).
6. Incubate at 37 °C, 5% CO₂ for **2-3 days**.

Day 5/6: Expansion

1. Centrifuge cells at 300 x g (RCF) for 5 min.
2. Aspirate supernatant.
3. Determine cell number.
4. Resuspend T cells at 1x10⁶ cells/ml in freshly prewarmed medium.
5. Add [recombinant human IL-2](#) (final concentration 20 IU/ml).
6. Incubate at 37 °C, 5% CO₂ for **2-3 days**.
7. Return to Day 1 for another cycle of activation/expansion.

This activation cycle can easily be performed up to 1 month with the same cell population, if the activation is repeated every 7-8 days.

RESTRICTION USE

This antibody is distributed for research purposes only. It is not intended for diagnosis or therapeutic use.

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

CULTURE CONDITIONS FOR T CELLS

Culture plate	Volume medium	Total cell number	bsAb CD3-CD28 (stock 0.1 mg/ml) to add per well
48 well	1 ml	1X10 ⁶	10 µl
24 well	2 ml	2X10 ⁶	20 µl
12 well	3 ml	3X10 ⁶	30 µl
6 well	5 ml	5X10 ⁶	50 µl
T25	10 ml	1X10 ⁷	100 µl
T75	20 ml	2X10 ⁷	200 µl

Table 1: Optimal surface density when working with purified T cells

RELATED PRODUCTS

Product	Description	Cat. Code
Recombinant human IL-2	Recombinant cytokine	rcyc-hil2
Normocin®	Antimicrobial reagent	ant-nr-1
Primocin®	Antimicrobial reagent	ant-pm-05

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