

**Description**

TL1A CHO Cell Line is a clonal CHO cell line expressing full-length human TL1A (TNF-like ligand 1A) (NM\_005118.4). The expressed TL1A may be released as a soluble form by proteolytic cleavage.

This cell line has been validated by flow cytometry, and in a co-culture assay with TL1A-Responsive Luciferase Reporter Jurkat Cell line (78811) in the presence of a TL1A neutralizing antibody.

**Background**

TL1A, also known as TNFSF15 (tumor necrosis factor superfamily member 15) or VEGI (Vascular endothelial growth inhibitor), is a member of the tumor necrosis factor family. It is expressed in different immune cells, such as monocytes, macrophages, dendritic cells, T cells, and non-immune cells. TL1A competitively binds to DR3 (death receptor 3), having a higher affinity for DcR3 (decoy receptor 3), and regulates proliferation, activation, apoptosis, and chemokine production in effector cells. The role of DR3 in T cell activation, and consequently in cytokine secretion and cell proliferation, makes it an attractive target in cancer therapy. Inhibition of DR3-TL1A interaction has substantial therapeutic potential in the treatment of solid tumors.

**Application(s)**

- Useful for cell-based TL1A binding assays.

**Materials Provided**

Components	Format
2 vials of frozen cells	Each vial contains $>1 \times 10^6$ cells in 1 ml of Cell Freezing Medium (BPS Bioscience #79796)

**Parental Cell Line**

CHO-K1 cells, Chinese Hamster Ovary, epithelial-like cells, adherent.

**Mycoplasma Testing**

The cell line has been screened to confirm the absence of Mycoplasma species.

**Materials Required but Not Supplied**

These materials are not supplied with the cell line but are necessary for cell culture and cellular assays. BPS Bioscience's reagents are validated and optimized for use with this cell line and are highly recommended for best results. Media components are provided in the Media Formulations section below.

**Media Required for Cell Culture**

Name	Ordering Information
Thaw Medium 3	BPS Bioscience #60186
Growth Medium 3J	BPS Bioscience #79974

*Materials Used in Cellular Assay*

Name	Ordering Information
Growth Medium 2A	BPS Bioscience #60190
Thaw Medium 2	BPS Bioscience #60184
TL1A Responsive Luciferase Reporter HEK293 Cell Line	BPS Bioscience #78811
Anti-TL1A Neutralizing Antibody	BPS Bioscience #101729
Tulisokibart	BPS Bioscience #83613
Clear-bottom, white 96-well tissue culture-treated plate	Corning #3610
ONE-Step™ Luciferase Assay System	BPS Bioscience #60690
Luminometer	

**Storage Conditions**

Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a -80°C freezer for long term storage. Contact technical support at [support@bpsbioscience.com](mailto:support@bpsbioscience.com) if the cells are not frozen in dry ice upon arrival.

**Media Formulations**

For best results, the use of validated and optimized media from BPS Bioscience is *highly recommended*. Other preparations or formulations of media may result in suboptimal performance.



Note: Thaw Media do *not* contain selective antibiotics. However, Growth Media *do* contain selective antibiotics, which are used to maintain selective pressure on the cell population expressing the gene of interest. Cells should be grown at 37°C with 5% CO<sub>2</sub>. BPS Bioscience's cell lines are stable for at least 10 passages when grown under proper conditions.

*Media Required for Cell Culture*

*Thaw Medium 3 (BPS Bioscience #60186):*

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin.

*Growth Medium 3J (BPS Bioscience #79974):*

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 5 µg/ml of Puromycin.

*Media required for Cellular Assay:*

*Thaw Medium 2 (BPS Bioscience #60184):*

RPMI1640 medium (ATCC modification) supplemented with 10% FBS, 1% Penicillin/Streptomycin.

*Growth Medium 2A (BPS Bioscience #60190):*

RPMI1640 medium (ATCC modification) supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 200 µg/ml Hygromycin B and 1 mg/ml G418.

## Cell Culture Protocol

### Cell Thawing

1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. As soon as the cells are thawed (it may be slightly faster or slower than 60 seconds), quickly transfer the entire contents of the vial to a tube containing 10 ml of pre-warmed Thaw Medium 3.

**Note: Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.**

2. Immediately spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed Thaw Medium 3.
3. Transfer the resuspended cells to a T25 flask or T75 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 48-72 hours of culture, check for cell attachment and viability. Change medium to fresh Thaw Medium 3 and continue growing in a 5% CO<sub>2</sub> incubator at 37°C until the cells are ready to passage.
5. Cells should be passaged before they reach 100% confluence. Switch to Growth Medium 3J for passage.

### Cell Passage

1. Aspirate the medium, wash the cells with phosphate buffered saline (PBS) without Ca<sup>2+</sup>/Mg<sup>2+</sup>, and detach the cells from the culture vessel with 0.25% Trypsin/EDTA, following the volumes recommended for the cell vessel being used.
2. Once the cells have detached, add Growth Medium 3J and transfer to a tube.
3. Spin down cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in Growth Medium 3J.
4. Seed into new culture vessels at the recommended sub-cultivation ratio of 1:6 to 1:8 once or twice a week.

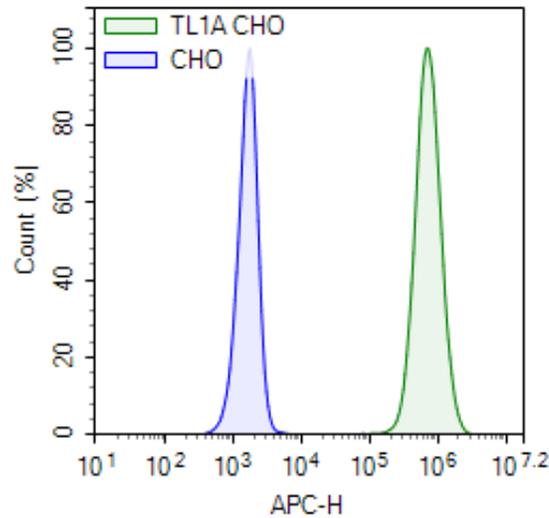
### Cell Freezing

1. After detachment, spin down the cells at 300 x g for 5 minutes.
2. Remove the medium and resuspend the cells in 4°C Cell Freezing Medium (BPS Bioscience #79796) at ~2 x 10<sup>6</sup> cells/ml.
3. Dispense 1 ml of cell suspension into each cryogenic vial. Place the vials in an insulated container for slow cooling and store at -80°C overnight.
4. Transfer the vials to liquid nitrogen the next day for long term storage.



Note: It is recommended to expand the cells and freeze at least 10 vials at an early passage for future use.

## Validation Data



*Figure 1: Cell surface expression of TL1A in TL1A CHO Cell Line assessed by flow cytometry.* TL1A CHO cells (green) and parental CHO-K1 (blue) were stained with the anti-TL1A antibody Tulisokibart (BPS Bioscience #83613) followed by Alexa Fluor® 647 anti-human IgG Fc Antibody (Jackson ImmunoResearch #109-606-008) and analyzed by flow cytometry. The y axis represents the cell %, while the x axis indicates Alexa Fluor 647 intensity.

## Functional Validation

### A. Inhibition of TL1A-induced reporter activity by Anti-TL1A Neutralizing Antibody, in a coculture assay with TL1A Responsive Luciferase Reporter Jurkat Cell Line.

- The following assay was designed for a 96-well format. To perform the assay in different tissue culture formats, the cell number and reagent volume should be scaled appropriately.
  - All conditions should be performed in triplicate.
  - We recommend using a non-specific antibody as control.
  - The assay should include “Antibody Treated Cells”, “Luminescence Background” and “Untreated Cells” conditions.
  - This assay requires TL1A Responsive Luciferase Reporter Jurkat Cell Line (#78811).
1. Thaw and grow TL1A Responsive Luciferase Reporter Jurkat cells in Thaw Medium 2 and Growth Medium 2A, respectively (for detailed information please refer to the datasheet of this cell line).

#### Day 1:

1. Prepare a cell suspension of TL1A CHO cells at  $2.0 \times 10^5$  cells/ml in Thaw Medium 3 (100  $\mu$ l/well).
2. Plate 100  $\mu$ l of TL1A CHO cells into each well of a white clear-bottom 96-well tissue culture plate. Leave a few wells empty (no cells) for the “Luminescence Background” control.

#### Day 2:

1. Prepare the test antibody at 2x the desired final concentrations in Thaw Medium 2 (50  $\mu$ l/well).

2. Prepare a cell suspension of TL1A Responsive Luciferase Reporter Jurkat cells in Thaw Medium 2 at  $8 \times 10^5$  cells/ml (50  $\mu$ l/well).
3. Remove the medium from TL1A CHO cells.
4. Add 50  $\mu$ l of diluted test antibodies to “Antibody Treated Cells” wells.
5. Add 50  $\mu$ l of Thaw Medium 2 to the “Untreated Cells” and “Luminescence Background” controls
6. Add 50  $\mu$ l of TL1A Responsive Luciferase Reporter Jurkat cells to the “Antibody Treated Cells”, and “Untreated Cells” wells.
7. Add 50  $\mu$ l of Thaw Medium 2 to the “Luminescence Background” control wells (for determining background luminescence).

*Note: If desired, TL1A CHO cells can be incubated with the anti-TL1A antibody for 30 minutes before TL1A Responsive Luciferase Reporter Jurkat cells are added to the wells for co-culture.*

8. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 5-6 hours.
9. Add 100  $\mu$ l of the ONE-Step™ Luciferase reagent per well.
10. Incubate with gentle agitation at RT for ~15 to 30 minutes.
11. Measure luminescence using a luminometer.
12. Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells. The fold induction of NF- $\kappa$ B luciferase reporter expression is the background-subtracted luminescence of “Antibody Treated Cells” well divided by the average background-subtracted luminescence of “Untreated Cells” control wells.

$$\text{Fold Induction} = \frac{\text{Lum}(\text{Antibody treated} - \text{Background})}{\text{Lum}(\text{Untreated} - \text{Background})}$$

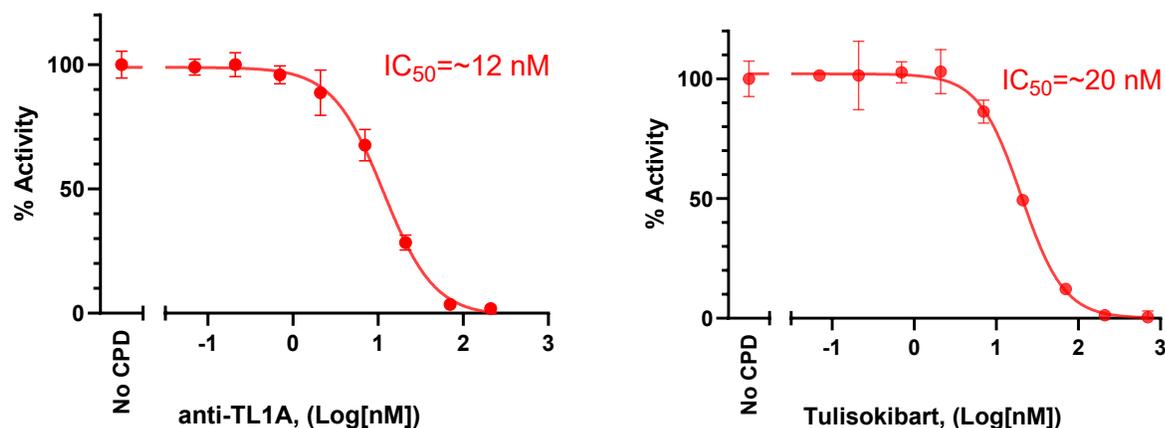


Figure 2: Inhibition of TL1A-induced reporter activity by Anti-TL1A Neutralizing Antibody, in the TL1A Responsive Luciferase Reporter Jurkat Cell Line cocultured with TL1A CHO Cell Line.

TL1A CHO cells were seeded overnight, and then co-cultured with TL1A Responsive Luciferase Reporter Jurkat Cells for 5 hours, in the presence of various amounts of Anti-TL1A Neutralizing Antibody (left, #101729) or Tulisokibart (right, #83613). TL1A-responsive luciferase activity was measured with ONE-Step™ Luciferase Assay System.

## References

Xu W. D., et al., 2022 *Front. Immunol.* 13: 1-10.  
Zwolak A., et al., 2022 *Sci. Rep.* 12(1): 20538.

## License Disclosure

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

## Troubleshooting Guide

Visit [bpsbioscience.com/cell-line-faq](https://bpsbioscience.com/cell-line-faq) for detailed troubleshooting instructions. For lot-specific information and all other questions, please visit <https://bpsbioscience.com/contact>.

## Related Products

Products	Catalog #	Size
Tulisokibart	83613	1 mg/ 5 mg
Anti-TL1A Neutralizing Antibody	101729	50 µg/ 100 µg
TL1A-Responsive Luciferase Reporter Jurkat Cell Line	78811	2 vials

Version 110325