



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

## **Data Sheet**

### ***CD70-CHO Recombinant Cell line***

**Cat. #: 79510**

#### **Product Description**

Recombinant CHO-K1 cells constitutively expressing human CD70 (also known as Tumor Necrosis Factor Ligand Superfamily Member 7, TNFSF7, CD27 Ligand, Ki-24 antigen, CD27-L, and CD27LG; GenBank accession #NM\_001252).

#### **Background**

CD70, also known as TNFSF7, is a cytokine which binds to CD27. CD70 is involved in T-cell activation and induces the proliferation of costimulated T-cells. Moreover, TNFSF7 enhances the generation of cytolytic T-cells. This cytokine is also reported to play a role in regulating B-cell activation, cytotoxic function of natural killer cells, and immunoglobulin synthesis. Diseases associated with CD70 include acute myocarditis, arthritis, and other inflammatory disorders.

#### **Applications**

This cell line is useful for CD70 binding assays, flow cytometry, or for screening for CD70 antibodies. Also useful for immune checkpoint co-stimulatory assay when co-cultured with CD27/NF- $\kappa$ B Reporter-Jurkat cell line (BPS Bioscience #79509).

#### **Format**

Each vial contains  $\sim 2 \times 10^6$  cells in 1 ml of 10% DMSO

#### **Storage**

Store in liquid nitrogen immediately upon receipt

#### **Mycoplasma Testing**

This cell line has been screened using the MycoAlert™ Mycoplasma Detection Kit (Lonza #LT07-118) to confirm the absence of Mycoplasma contamination.

#### **Culture Medium**

**Thaw Medium 3 (BPS Cat #60186):** F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin.

**Growth Medium 3D (BPS Bioscience, #79539):** F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 1 mg/ml G418.

Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 3D.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

### Recommended Culture Condition

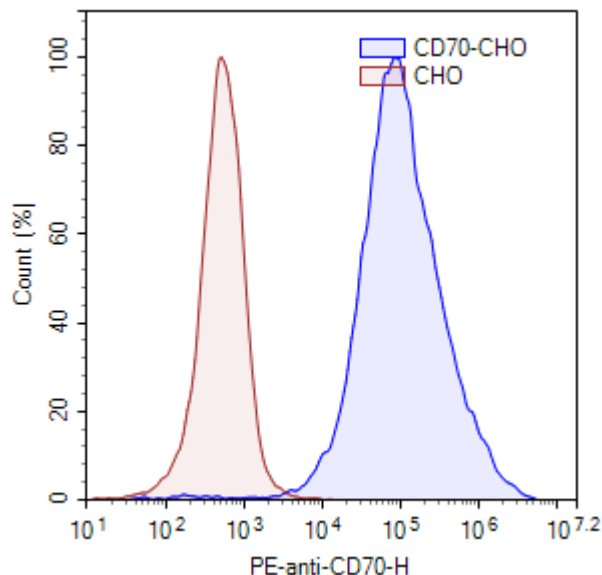
It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 3 (**no G418**), spin down cells, re-suspend cells in pre-warmed Thaw Medium 3 (**no G418**), transfer re-suspended cells to T25 flask and culture in 37°C CO<sub>2</sub> incubator overnight. The next day, replace the medium with fresh Thaw Medium 3 (**no G418**), and continue growing culture in a CO<sub>2</sub> incubator at 37°C until the cells are ready to be split. Cells should be split before they reach complete confluence. At first passage switch to Growth Medium 3D (**contains 1 mg/ml G418**).

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add Growth Medium 3D and transfer to a tube, spin down cells, re-suspend cells and seed appropriate aliquots of cell suspension into new culture vessels.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 0.05% Trypsin/EDTA. After detachment, add Thaw Medium 3 (**no G418**) and count the cells, then transfer to a tube, spin down cells, and resuspend in Freezing Medium (10% DMSO + 90% FBS) at  $\sim 2 \times 10^6$  cells/ml. Dispense 1 ml of cell aliquots into cryogenic vials. Place vials in an insulated container for slow cooling and store at -80°C overnight. Transfer to liquid nitrogen the next day for storage. It is recommended to expand the cells and freeze down more than 10 vials of cells for future use at early passage.

### Functional Validation and Assay Performance

Expression of human CD70 in CHO-K1 cells was confirmed by FACS.



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

## Sequence

Human CD70 sequence (accession number NM\_001252)

MPEEGSGCSVRRRPYGCVLRAALVPLVAGLVICLVVCIQRFAQAQQQLPLESLGWDVAELQLNHT  
GPQQDPRLYWQGGPALGRSFLHGPELDKGQLRIHRDGIYMVHIQVTLAICSSTTASRHHPTTLAV  
GICSPASRSISLLRLSFHQGCTIASQRLTPLARGDTLCTNLTGTLTLLPSRNTDETFFGVQWVRP

## Related Products

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
CD27, Fc fusion Protein	71176	100 µg
CD70(CD27L), His-tag Protein	71178	100 µg
CD27 CHO-K1 Stable Recombinant Cell Line	60624	2 vials
CD27/NF-κB Reporter-Jurkat Recombinant Cell Line	79510	2 vials
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
Thaw Medium 2	60184	100 ml
Thaw Medium 3	60186	100 ml

### License Disclosure:

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

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)