



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### **TCR Activator - CHO Recombinant Cell line Cat#: 60539**

#### **Product Description**

Recombinant CHO-K1 cell constitutively expressing a membrane bound, engineered T cell receptor (TCR) activator.

#### **Application**

- *in vitro* stimulation of T cells.
- Control cell line for TCR activator / PD-L1 - CHO cell line, BPS Cat. #60536

#### **Format**

Each vial contains 2.5 X 10<sup>6</sup> cells in 1 ml of 10% DMSO

#### **Storage**

Immediately upon receipt, store in liquid nitrogen.

#### **Mycoplasma Testing**

The cell line has been screened using the PCR-based Venor<sup>®</sup>GeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of *Mycoplasma* species.

#### **General Culture Conditions**

**Thaw Medium 3 (BPS Cat. #60186):** Ham's F-12 medium (Hyclone # SH30526.01) supplemented with 10% FBS (Life technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

**Growth Medium 3B (BPS Cat. #79529):** Thaw Medium 3 plus 500 µg/ml of Hygromycin B (Hyclone #SV30070.01) to ensure the recombinant expression.

Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 3B. TCR activator - CHO cells should exhibit a typical cell division time of ~24 hours.

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 Hygromycin B**), spin down the cells, and resuspend the cells in pre-warmed Thaw Medium 3 (**no Hygromycin B**). Transfer resuspended cells to a T-25 flask and culture at 37°C in a CO<sub>2</sub> incubator overnight. The next day, replace the medium with fresh Thaw Medium 3 (**no Hygromycin B**), and continue growing culture in a CO<sub>2</sub> incubator at 37°C until the cells

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

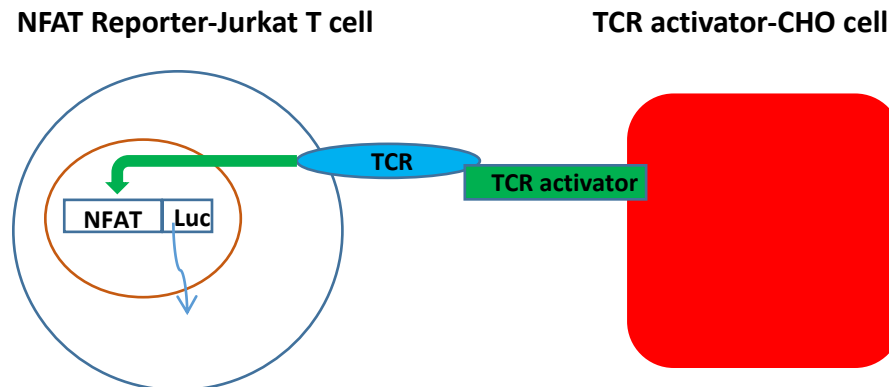
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

are ready to be split. Cells should be split before they reach complete confluence. At first passage, switch to Growth Medium 3B (**contains Hygromycin B**).

To passage the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 0.05% Trypsin/EDTA. Add Growth Medium 3B and transfer to a tube, and spin down the cells. Re-suspend the cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ration: 1:10 to 1:20 twice a week.

### Functional Validation and Assay Performance

The functionality of the cell line was validated using a luciferase reporter cell-based assay. In this assay, Jurkat T cells expressing NFAT reporter are co-cultivated with TCR activator – CHO cells. TCR complexes on Jurkat cells are activated by TCR activator on TCR activator – CHO cells, resulting in expression of the NFAT luciferase reporter.



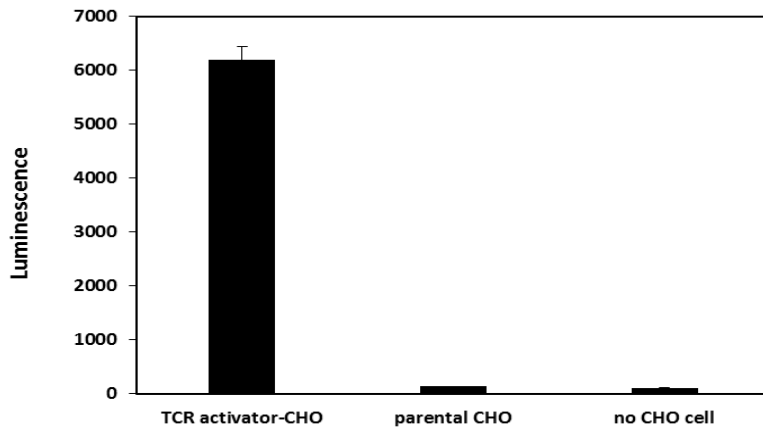
**Figure 1. Co-culture of TCR activator-CHO cells with NFAT Reporter-Jurkat cells induced NFAT reporter expression in Jurkat cells.**

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

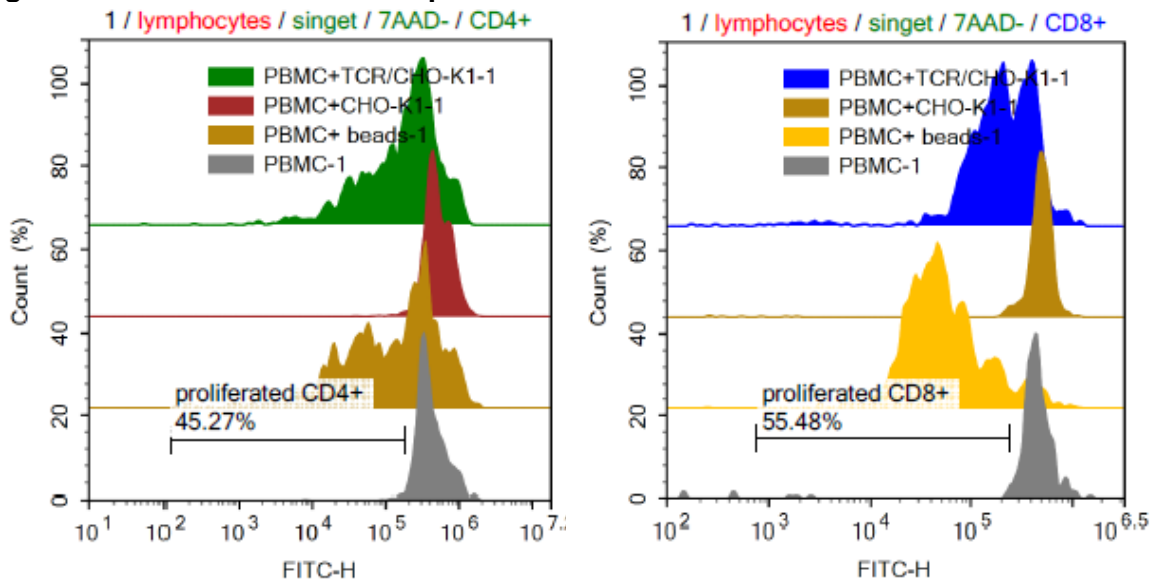
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



TCR activator-CHO cells or parental CHO cells were seeded in 96-well plate. The next day, NFAT Reporter-Jurkat cells (BPS Cat. #60621) were added to TCR activator-CHO cells. After incubation, ONE-Step™ Luciferase reagent (BPS Cat. #60690) was added to the cells to measure NFAT activity.

**Figure 2. TCR activator/CHO-K1 promotes proliferation of T cells.**



PBMCs were stained with CellTrace™ CFSE (Thermo Fisher Cat # C34554) and co-cultured with TCR-activator CHO-K1, wildtype CHO-K1 (PBMC+CHO-K1), activated with anti-CD3/CD28 beads (PBMC + beads), or untreated (PBMC) for 72 hours. (Left) CD4+ proliferation ; (Right) CD8+ proliferation.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Related Products

<b><u>Product</u></b>	<b><u>Cat. #</u></b>
NFAT Reporter – Jurkat cell line	60621
PD-1/NFAT Reporter-Jurkat cell line	60535
TIGIT/NFAT Reporter – Jurkat cell line	60538
TCR activator / PD-L1 - CHO cell line	60536
TCR activator / PD-L1 expression kit	60610
TCR activator / PD-L2 expression kit	60620
ONE-Step™ Luciferase Assay System	60690

## **License Disclosure:**

Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use, and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact [sales@bpsbioscience.com](mailto:sales@bpsbioscience.com) for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

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