



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

TCR Activator - Raji Cell Line

Catalog #: 60556

Product Description

Recombinant Raji cell expressing an engineered T cell receptor (TCR) activator.

Format

Each vial contains 1×10^6 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[®]GeM Mycoplasma Detection kit (Sigma-Aldrich, #MP0025) to confirm the absence of *Mycoplasma* species.

Culture Medium and Recommended Culture Conditions

Thaw Medium 2 (BPS Cat. #60184): RPMI1640 medium (Life Technologies, #A10491-01) supplemented with 10% FBS (Life Technologies, #26140-079), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01).

Complete Growth Medium: Thaw Medium 2 (BPS Bioscience, #60184) + 200 µg/ml of Hygromycin B (Hyclone, #SV30070.01).

Cells should be grown at 37°C with 5% CO₂ using **complete growth medium** (Thaw Medium 2 + 200 µg/ml of Hygromycin B).

To thaw the cells, it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath. Transfer the entire contents of the vial to a tube containing 10 ml of Thaw Medium 2 (**no Hygromycin B**), spin down cells at 1500 rpm, remove supernatant and resuspend cells in 5 ml of pre-warmed Thaw Medium 2 (**no Hygromycin B**). Transfer resuspended cells to a T25 flask and culture at 37°C in a 5% CO₂ incubator overnight. The next day, add an additional ~3 ml of fresh warm Thaw Medium 2 (**no Hygromycin B**), and continue growing culture in a CO₂ incubator at 37°C until the cells are ready to be split. Cells should be split before they reach 2.5×10^6 cells/ml. At first passage, switch to complete growth medium (**contains Hygromycin B**).

To passage the cells, dilute cell suspension into new culture vessels at no less than 0.2×10^6 cells/ml. Subcultivation ratio: ~1:5 to 1:10 once or twice a week, so cells are maintained at 0.2×10^6 cells/ml to 2.5×10^6 cells/ml.

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

Please visit our website at: www.bpsbioscience.com

***Note:** Just after thawing, the cells may grow at a slower rate. It is recommended to split the cells at no less than 0.4×10^6 cells/ml at the beginning of culturing. After two or three passages, the cell growth rate increases and the cells can be split to 0.2×10^6 cells/ml.*

To freeze down the cells, spin down cells, and resuspend cell pellet in 4°C Freezing Medium (10% DMSO + 90% FBS) to $\sim 2 \times 10^6$ 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 at early passage, freeze down more than 10 vials of cells for future use.

Functional Validation

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 – Raji cells. TCR complexes on the Jurkat cells are activated by TCR activator on the TCR Activator – Raji cells, resulting in expression of the NFAT luciferase reporter.

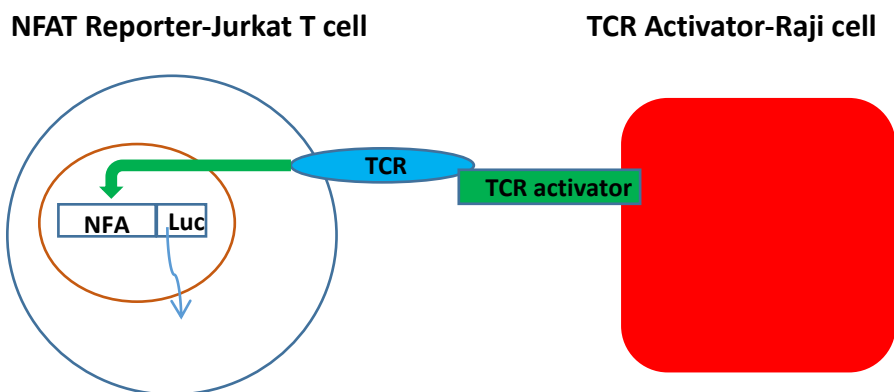


Figure 1. Co-culture of TCR Activator-Raji cells with NFAT Reporter-Jurkat cells induced NFAT reporter expression in Jurkat cells.

TCR Activator-Raji cells or parental Raji cells were seeded in 96-well plate, then NFAT Reporter-Jurkat cells (BPS Bioscience, #60621) were added to TCR activator-Raji cells. After 5-6 hour incubation, ONE-Step™ Luciferase reagent (BPS Bioscience, #60690) was added to the cells to measure NFAT-induced luciferase activity.

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**

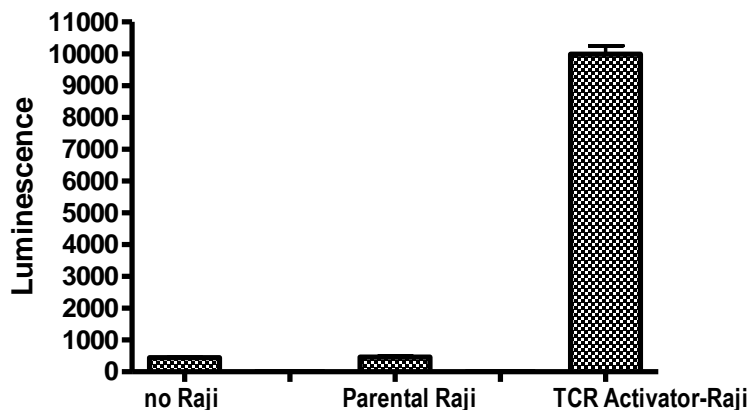
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: 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

NFAT Reporter-Jurkat



Product

TCR Activator-CHO cell line
NFAT Reporter – Jurkat cell line

Thaw Medium 2

ONE-Step™ Luciferase Assay System

ONE-Step™ Luciferase Assay System

TCR Activator / PD-L1 Mammalian Expression Kit

TCR Activator / PD-L2 Mammalian Expression Kit

PD-L1 / TCR Activator - CHO Recombinant Cell Line

CD155 / TCR Activator - CHO Recombinant Cell line

PD-1 / NFAT Reporter - Jurkat Recombinant Cell Line

TIGIT / NFAT Reporter - Jurkat Recombinant Cell Line

LAG3 / NFAT Reporter - Jurkat Recombinant Cell Line

Cat. #

60539

60621

60184

60690-1

60690-2

60610

60620

60536

60548

60535

60538

71278

Size

2 vials

2 vials

100 ml

10 ml

100 ml

500 rxns.

500 rxns.

2 vials

2 vials

2 vials

2 vials

2 vials

Notes

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 the 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 Bioscience 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 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

Please visit our website at: www.bpsbioscience.com