



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### **ICOS/NFAT Reporter-Jurkat Recombinant Cell Line**

### **Catalog #: 79668**

#### **Product Description**

Recombinant Jurkat T cell expressing firefly luciferase gene under the control of NFAT response elements with constitutive expression of a chimeric receptor consisting human ICOS (also known as inducible T-cell costimulator or CD278, Genbank Accession #NM\_012092) and the cytoplasmic domain of human CD3 zeta.

#### **Background**

ICOS is a costimulatory molecule of the CD28 cell surface receptor superfamily that is expressed on activated T-cells. ICOS is involved in T-cell responses upon binding with its ligand, ICOSL (also known as B7-H2, CD275), which is normally expressed on B-cells, dendritic cells and monocytes. ICOS expression confers an activated phenotype and a strong suppressive capacity to intra-tumoral regulatory T-cells. The ICOS/ICOSL pathway is a key target for cancer immunotherapy.

#### **Application**

- Screen for agonists of ICOS signaling in a physiologically relevant cellular context
- Characterize T cell-mediated immune responses of ICOS and its interaction with ICOSL

#### **Format**

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

#### **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.

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

**Growth Medium 2F (BPS Bioscience #79669):** Thaw Medium 2 (BPS Bioscience #60184), 1 mg/ml of Geneticin (Life Technologies #11811031), and 0.5 µg/ml of Puromycin (Takara, #631306).

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

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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

### Recommended Culture Condition

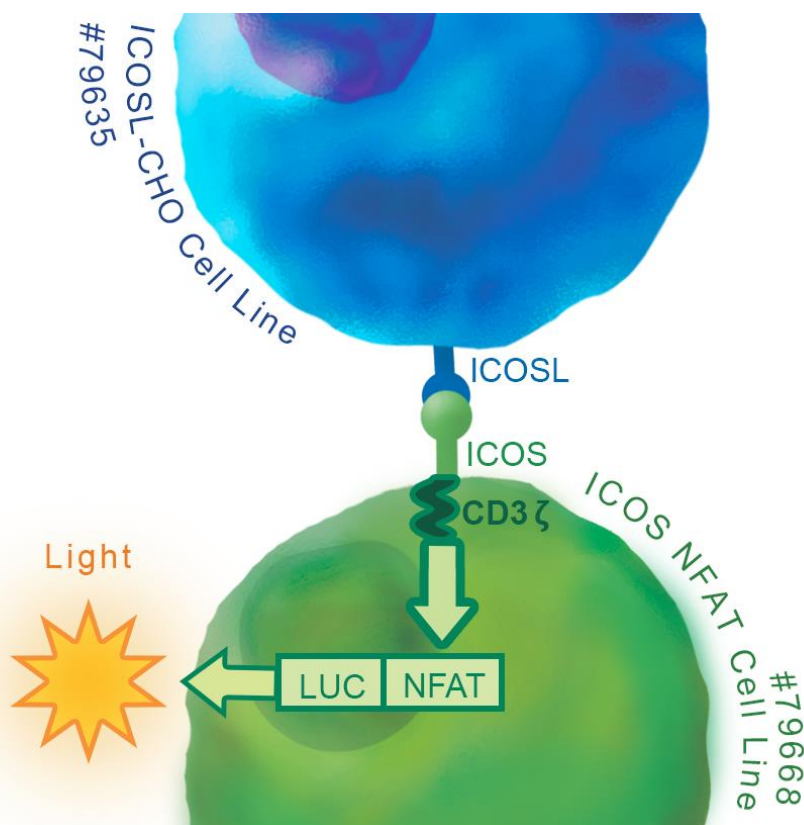
It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, then transfer the entire contents of the vial to a tube containing 10 ml of Thaw Medium 2 (no Geneticin and Puromycin). Spin down the cells, remove supernatant and resuspend cells in pre-warmed Thaw Medium 2 (no Geneticin and Puromycin). Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator. After 24 hours of culture, add an additional 3 – 4 ml of growth medium without antibiotics. At first passage, switch to growth medium 2F (contains Geneticin and Puromycin). Cells should be split before they reach 2 x 10<sup>6</sup> cells/ml. Note: This cell line tends to grow more slowly than the parental Jurkat cells.

To passage the cells, dilute cell suspension into new culture vessels at no less than 0.2 x 10<sup>6</sup> cells/ml. Subcultivation ratio: 1:10 to 1:20 twice a week.

### Functional Validation and Assay Performance

Expression of human ICOS in the Jurkat cell line was confirmed by Flow Cytometry. The functionality of the cell line was validated using an anti-ICOS agonist antibody and a ICOS:ICOSL cell-based assay.

### Assay Principle



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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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

### Materials Required but Not Supplied

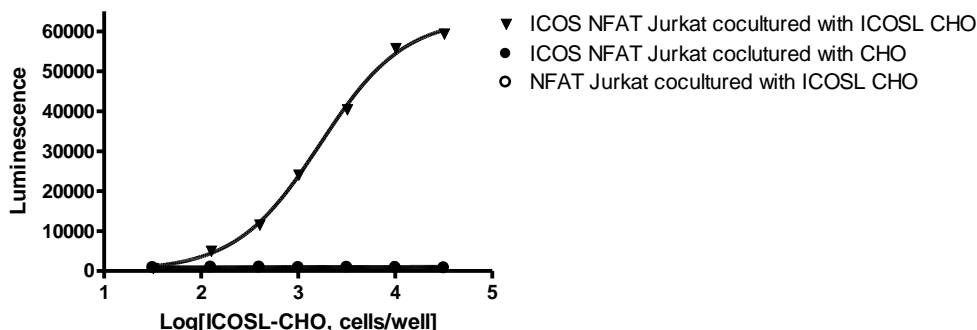
- Assay Medium: Thaw Medium 2 (BPS Bioscience #60184)
- ICOSL-CHO K1 Recombinant Cell Line (BPS Bioscience #79635) and its recommended growth medium (Growth Medium 3D, BPS Bioscience #79539)
- 96-well tissue culture-treated white clear-bottom assay plate
- One-Step luciferase assay system (BPS Bioscience #60690) for measuring firefly luciferase activity
- Luminometer

### Assay Protocol and Functional Analysis

#### A) ICOS/NFAT reporter activities stimulated by ICOSL-CHO cells

1. Harvest ICOSL-CHO cells from culture in growth medium and seed cells into white clear-bottom 96-well microplate in 100  $\mu$ l growth medium. Incubate the plate at 37°C in a CO<sub>2</sub> incubator overnight. Leave a few wells empty as cell-free controls.
2. Next day, harvest the ICOS/NFAT Reporter-Jurkat cells by centrifugation and resuspend in assay medium. Dilute the cells to 3 x 10<sup>5</sup> / ml in assay medium. Remove the medium from ICOSL-CHO cells and add 50  $\mu$ l of ICOS/NFAT Reporter-Jurkat cells to the wells.
3. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 5 hours.
4. Perform luciferase assay using the ONE-Step luciferase assay system: Prepare the ONE-Step Luciferase reagent per recommended protocol. Add 50  $\mu$ l of ONE-Step Luciferase reagent per well and rock at room temperature for ~30 minutes. Measure luminescence using a luminometer.
5. Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells.

**Figure 1. The ICOS/NFAT Reporter Activities Stimulated by ICOSL-CHO Cells.**



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

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

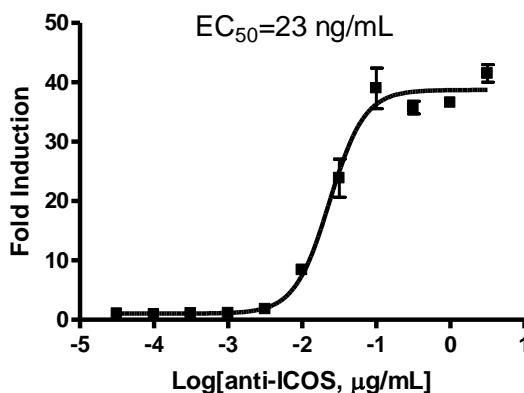
**B) Functional assay of anti-ICOS agonist antibody on ICOS/NFAT Reporter-Jurkat cells**

1. Harvest ICOS/NFAT Reporter-Jurkat cells from culture in growth medium and seed 20,000 cells per well into white clear-bottom 96-well plate in 45  $\mu$ l of assay medium.
2. Dilute anti-ICOS agonist antibody in assay medium and treat cells with 5  $\mu$ l of 10X dilutions of the anti-ICOS antibody.
3. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for 5 hours.
4. Perform luciferase assay using the ONE-Step luciferase assay system: Prepare the ONE-Step Luciferase reagent per recommended protocol. Add 50  $\mu$ l of ONE-Step Luciferase reagent per well (BPS Bioscience #60690) and rock at room temperature for 20 minutes. Measure luminescence using a luminometer

Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells. The fold induction of NFAT luciferase reporter expression = background-subtracted luminescence of stimulated well / average background-subtracted luminescence of unstimulated control wells.

**Figure 2 Dose response of anti-ICOS agonist antibody on ICOS/NFAT Reporter-Jurkat cells**

Serial dilutions of anti-ICOS agonist antibody (Biolegend #313512) were added to ICOS/NFAT Reporter-Jurkat cells (BPS Bioscience #79668), and then incubated at 37°C for 5 hours. After the treatment, the Luciferase assay was performed.



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

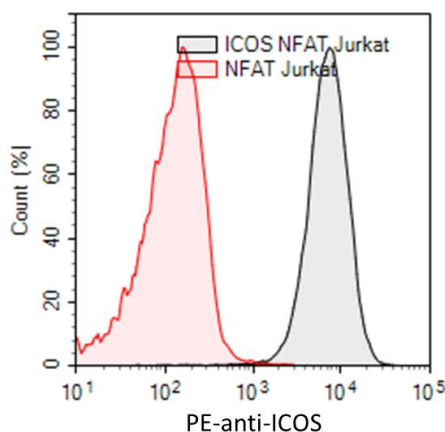
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

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

**Figure 3. FACS Analysis of Cell Surface Expression of ICOS in ICOS/NFAT Reporter-Jurkat cells.**

ICOS/NFAT Reporter-Jurkat cells or NFAT Reporter-Jurkat cells (BPS Bioscience #60621) were stained with PE-labeled anti-ICOS antibody (BD Bioscience #557802, clone#DX29) and analyzed by FACS. Y-axis is the cell count. X-axis is the intensity of PE.



**License Disclosure:**

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

**Related Products**

<b>Related Products</b>	<b>Cat. #</b>	<b>Size</b>
ICOSL (B7-H2) -CHO Recombinant Cell Line	79635	2 vials
NFAT Reporter (Luc) – Jurkat Recombinant Cell Line	60621	2 vials
ICOS (CD278), Fc fusion (Human)	71179	100 µg
B7-H2 (CD275, ICOSL), Fc fusion (Human) HiP™	71130	100 µg
B7-H2, Avi-His-Tag	79119	100 µg
B7-H2, Avi-His-Tag, Biotin-Labeled	79300	50 µg
CD276 (B7-H3), Avi-His-Tag HiP™	79337	100 µg
B7-H4, His-tag (Human)	71144	100 µg
Thaw Medium 2	60184	100 ml
Growth Medium 2F	79669	500 ml
Growth Medium 3D	79539	500 ml

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

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