



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

STAT5 Reporter (Luc) – Ba/F3 Cell line

Catalog #: 79772

Product Description

The STAT5 Reporter (Luc)-Ba/F3 cell line is designed for monitoring STAT5 signal transduction pathways. It contains a firefly luciferase gene driven by the STAT5 response element located upstream of the minimal TATA promoter. After activation by cytokines or growth factors, endogenous STAT5 binds to the DNA response elements, inducing transcription of the luciferase reporter gene.

Application

- Monitor STAT5 signaling pathway activity
- Screen for activators or inhibitors of STAT5 signaling pathway

Format

Each vial contains ~3 x 10⁶ cells in 1 ml of FBS with 10% DMSO

Storage

Immediately upon receipt, store in liquid nitrogen.

Host Cell

Ba/F3: IL-3 dependent murine pro B cell line. Non-adherent cells.

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.

General Culture Conditions

Thaw Medium 8 (BPS Bioscience, #79652) plus 5 ng/mL mouse IL-3: RPMI1640 medium (Life Technologies, #A10491-01) supplemented with 10% heat-inactivated FBS (Life Technologies, #10082147), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01), 5 ng/ml mIL-3 (R&D systems, #403-ML-010).

Growth Medium 8A (BPS Bioscience, #79653) plus 5 ng/mL mouse IL-3: Thaw Medium 8 (BPS Bioscience, #79652), 1 µg/ml of Puromycin (Takara, #631306), 5 ng/ml mIL-3 (R&D systems, #403-ML-010).

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 8A with the addition of 5 ng/ml mIL-3. **Note:** mIL-3 is essential for Ba/F3 cell maintenance. Thaw Medium 8 and Growth Medium 8A do not contain mIL-3.

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

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

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 8 plus 5 ng/ml mIL-3 (**no Puromycin**). Spin down the cells, remove supernatant and resuspend cells in pre-warmed Thaw Medium 8 plus 5 ng/ml mIL-3 (**no Puromycin**). Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO₂ incubator. After 24 hours of culture, add an additional 1 – 2 ml of Thaw Medium 8 plus 5 ng/ml mIL-3 (**no Puromycin**). At first passage, switch to Growth Medium 8A (**contains Puromycin**) plus 5 ng/ml IL-3. Cells should be split before they reach 2.0 x 10⁶ cells/ml.

To passage the cells, dilute cell suspension into new culture vessels at no less than 0.1 x 10⁶ cells/ml. Do not allow the cell density to exceed 2.0 x 10⁶ cells/ml.

Assay Performance

The following assays are designed for 96-well format. To perform the assay in different tissue culture formats, cell number and reagent volume should be scaled appropriately.

Materials Required but Not Supplied

- mIL-3 (R&D systems, #403-ML-010)
- Assay Medium: Thaw Medium 8 (BPS Bioscience, #79652)
- Growth Medium 8A (BPS Bioscience, #79653)
- 96-well tissue culture treated white clear-bottom assay plate (Corning, #3610)
- One-Step luciferase assay system (BPS Bioscience, #60690)
- Luminometer

mIL-3 Dose Response

1. Harvest STAT5 Reporter (Luc) Ba/F3 cells from culture in Growth Medium 8A containing 5 ng/ml mIL-3. Wash cells once with assay medium (containing no mIL-3); resuspend the cells in assay medium and seed cells into a white opaque 96-well microplate at a density of ~20,000 cells per well in 90 µl of assay medium. Leave a couple wells empty for use as a background control.
2. Prepare threefold serial dilution of mIL-3 in assay medium. Add 10 µl of diluted mIL-3 to mIL-3 stimulated wells.
3. Add 10 µl of assay medium to the unstimulated control wells (for measuring the uninduced level of STAT5 reporter activity).
4. Add 100 µl of assay medium to the cell-free control wells (for determining background luminescence).
5. Incubate at 37°C with 5% CO₂ for 5-16 hours.
6. Prepare ONE-Step™ Luciferase Assay reagent per recommended instructions. Add 100 µl of ONE-Step™ Luciferase reagent per well. Incubate at room temperature for ~15 to 30

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**

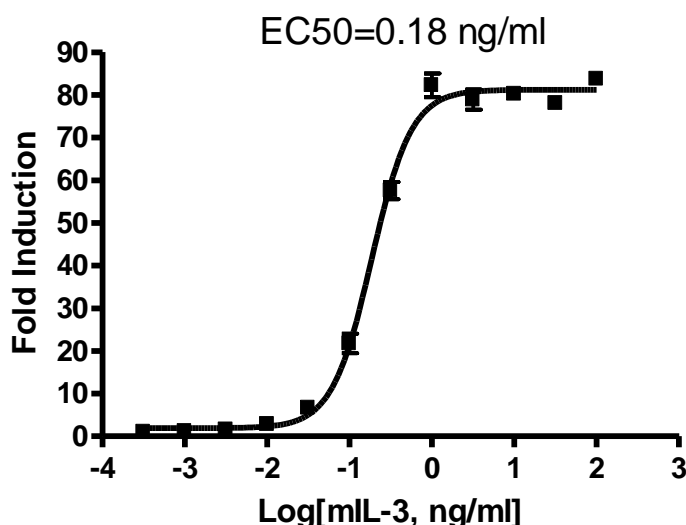
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

minutes and measure luminescence using a luminometer. Subtract background luminescence value from all readings.

Figure 1. mL-3 dose response in STAT5 reporter (Luc)-Ba/F3 cells. Cells were treated with mL-3 for ~ 16 hours. The results were shown as fold induction of luciferase reporter expression. Fold induction was determined by comparing values against the mean value for control cells without mL-3 treatment.

The EC50 of mL-3 in this cell line is ~0.18 ng/ml.



Related Products

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
Thaw Medium 8	79652	100 ml
Growth Medium 8A	79653	500 ml
Mouse IL-3	403-ML-010	2 µg
Mouse IL-3	403-ML-010	10 µg

References

1. Palacios, R., Henson, G., Steinmetz, M., McKearn, J. P. (1984). Interleukin-3 supports growth of mouse pre-B-cell clones in vitro. *Nature*, **309 (5964)**: 126-131.
2. Jaster R, Tschirch E, Bittorf T, Brock J. (1999). Role of STAT5 in interferon-alpha signal transduction in Ba/F3 cells. *Cell Signal.*, **11(5)**:331-5.

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

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

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 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com