



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

NF-κB Reporter (Luc) – NIH/3T3 Cell Line

Catalog #: 79469

Product description

The NF-κB reporter (Luc)-NIH/3T3 cell line is designed for monitoring nuclear factor Kappa B (NF-κB) signal transduction pathways. It contains a firefly luciferase gene driven by four copies of the NF-κB response element located upstream of the minimal TATA promoter. After activation by pro-inflammatory cytokines or stimulants of lymphokine receptors, endogenous NF-κB transcription factors bind to the DNA response elements, inducing transcription of the luciferase reporter gene.

Application

- Monitor NF-κB signaling pathway activity.
- Screen for activators or inhibitors of NF-κB signaling pathway.

Format

Each vial contains ~2 x 10⁶ 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.

General Culture Conditions

Thaw Medium 5 (BPS Cat. #60182): DMEM (Hyclone #SH30243.01), supplemented with 10% Calf Bovine Serum (Hyclone #SH3007203), 1% Penicillin/Streptomycin (Hyclone #SV30010.01).

Growth Medium 5A (BPS Cat. #79534): Thaw Medium 5 (BPS Cat. #60186) and 600 µg/ml of Geneticin (Life Technologies #11811031).

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 5A.

NF-κB reporter (Luc) – NIH/3T3 cells should exhibit a typical cell division time of 24 hours.

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

To thaw the cells, 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 5 (**no Geneticin**). Spin down the cells, remove supernatant and resuspend cells in pre-warmed Thaw Medium 5 (**no Geneticin**). Transfer the resuspended cells to a T75 flask and incubate at 37°C in a 5% CO₂ incubator. After 24 hours of culture, add an additional 3 – 4 ml of Thaw Medium 5 (**no Geneticin**). At first passage, switch to Growth Medium 5A (**contains Geneticin**).

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add Growth Medium 5A and transfer to a tube. Spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:5 to 1:10, twice a week.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with Trypsin/EDTA. Add Growth Medium 5A and transfer to a tube, spin down cells, and resuspend in freezing medium (10% DMSO + 90% FBS). Place at -80°C overnight and place in liquid nitrogen the next day. Alternatively, vials may be placed directly in liquid nitrogen.

Assay Performance

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

Materials Required but Not Supplied

- Mouse TNF α (BioLegend #754402-10 μ G)
- Assay medium: Thaw Medium 5 (BPS Bioscience #60182) or
 - DMEM medium (Hyclone #SH30243.01) + 10% Calf Bovine Serum (Hyclone #SH3007203) + 1% Pen/Strep (Hyclone #SV30010.01)
- Growth Medium 5A (BPS Cat. #79534)
- 96-well tissue culture treated white clear-bottom assay plate (Corning #3610)
- One-Step luciferase assay system (BPS Bioscience #60690)
- Luminometer

A. TNF α dose response

1. Harvest NF- κ B reporter (Luc)-NIH/3T3 cells from culture in Growth Medium 5A and seed cells at a density of ~25,000 cells per well into white opaque 96-well microplate in 50 μ l of assay medium. Incubate cells at 37°C with 5% CO₂ overnight.
2. Prepare threefold serial dilution of TNF α in assay medium. Add 50 μ l of diluted TNF α to TNF α -stimulated wells.

Add 50 μ l of assay medium to the unstimulated control wells (for measuring uninduced level of NF- κ B reporter 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

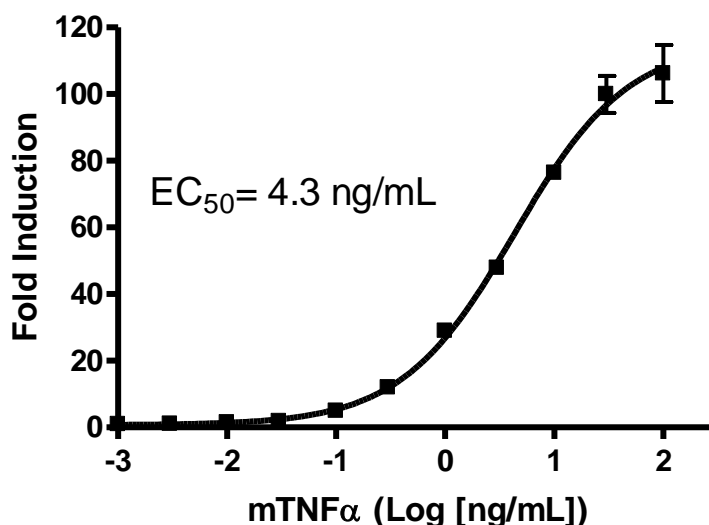
Please visit our website at: www.bpsbioscience.com

Add 100 μ l of assay medium to cell-free control wells (for determining background luminescence).

5. Incubate at 37°C with 5% CO₂ for ~3-6 hours.
6. Prepare ONE-Step™ Luciferase Assay reagent as directed. Add 100 μ l of ONE-Step™ reagent per well. Incubate at room temperature for ~15 to 30 minutes and measure luminescence using a luminometer.
7. Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells. The fold induction of NF- κ B luciferase reporter expression = background-subtracted luminescence of stimulated well / average background-subtracted luminescence of unstimulated control wells.

Figure 1. TNF α dose response in NF- κ B reporter (Luc)-NIH/3T3 cells. Cells were treated with TNF α for ~ 6 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 TNF α treatment.

The EC₅₀ of mTNF α in this cell line is ~4.3 ng/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



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

Related Products

| Product | Cat. # | Size |
|---|---------------|-------------|
| NF-κB reporter (Luc) - HEK293 Cell line | 60650 | 2 vials |
| NF-κB reporter (Luc) - Jurkat Cell line | 60651 | 2 vials |
| NF-κB reporter (Luc) - CHOK1 Cell line | 60622 | 2 vials |
| NF-κB reporter (Luc) – A549 Cell line | 60625 | 2 vials |
| ONE-Step™ Luciferase Assay System | 60690-1 | 10 ml |
| ONE-Step™ Luciferase Assay System | 60690-2 | 100 ml |
| NF-κB Reporter Kit (NF-κB Signaling Pathway) | 60614 | 500 rxns. |
| Transfection Collection™ : | | |
| NF-κB Transient Pack (NF-κB Signaling Pathway) | 79268 | 500 rxns. |
| Transfection Collection™ : | | |
| NF-κB Reporter Cellular Assay Pack (HEK293) | 79327 | 2 vials |
| Transfection Collection™ : | | |
| NF-κB Reporter Cellular Assay Pack (HCT116) | 79326 | 2 vials |

References

1. Pessara U, Koch N (1990) Tumor necrosis factor alpha regulates expression of the major histocompatibility complex class II-associated invariant chain by binding of an NF-κB-like factor to a promoter element. *Mol Cell Biol.* **10(8)**:4146-4154.
2. Baeuerle PA (1998) Pro-inflammatory signaling: last pieces in the NF-κB puzzle? *Curr Biol.* **8(1)**:R19-R22.

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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com