



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- κ B Reporter (Luc) - Jurkat Cell line **Catalog #: 60651**

Product Description

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

Application

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

Format

Each vial contains $\sim 2 \times 10^6$ 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) to confirm the absence of *Mycoplasma* species.

General Culture Conditions

Thaw Medium 2 (BPS Bioscience #60184): RPMI 1640 medium (Thermo Fisher, Cat. #A1049101) supplemented with 10% FBS (Thermo Fisher, Cat. #26140079), 1% Penicillin/Streptomycin (Hyclone #SV30010.01).

Growth Medium 2B (BPS Bioscience #79530): Thaw Medium 2 plus 1 mg/ml of Geneticin (Thermo Fisher, Cat. #11811031).

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 2B.

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**). Spin down the cells, remove supernatant and resuspend cells in pre-warmed Thaw Medium 2, (**no Geneticin**). 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 3 – 4 ml of Thaw Medium 2, (**no Geneticin**). At first passage, switch to Growth Medium 2B (**contains geneticin**). Cells should be split before they reach 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



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 passage the cells, dilute cell suspension into new culture vessels at no less than 0.1×10^6 cells/ml. Subcultivation ratio: 1:5 to 1:10 twice a week.

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

- TNF α (Sigma, #T0157-10UG)
- PMA (LC Laboratories, #P1680)
- Ionomycin (Sigma, #I3909)
- Assay Medium: Thaw Medium 2 (BPS Bioscience, #60184)
- Growth Medium 2B (BPS Bioscience, #79530)
- 96-well tissue culture treated white clear-bottom assay plate (Corning # 3610)
- One-Step luciferase assay system (BPS Bioscience, #60690) or other luciferase reagents for measuring firefly luciferase activity
- Luminometer

A. TNF α dose response

1. Harvest NF- κ B reporter (Luc)-Jurkat cells from culture in Growth Medium 2B and seed cells at a density of $\sim 40,000$ cells per well into white clear-bottom 96-well assay plate 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.
3. Add 50 μ l of assay medium to the unstimulated control wells (for measuring uninduced level of NF- κ B reporter activity).
4. 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. Add 100 μ l of ONE-Step™ Luciferase Assay reagent per well. Incubate at room temperature for ~ 15 to 30 minutes and measure luminescence using a luminometer.

Figure 1. TNF α dose response in NF- κ B reporter (Luc)-Jurkat cells. Cells were treated with TNF α for ~ 6 hours. The results were shown as fold induction of luciferase reporter expression.

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

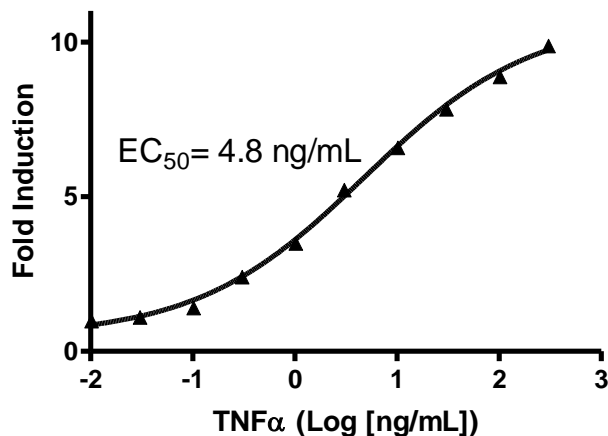
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

Fold induction was determined by comparing values against the mean value for control cells without TNF α treatment.

The EC₅₀ of TNF α in this cell line is ~4.8 ng/ml.



B. PMA dose response

1. Harvest NF- κ B reporter (Luc)-Jurkat cells from culture in Growth Medium 2B and seed cells at a density of ~40,000 cells per well into white clear-bottom 96-well assay plate in 50 μ l of assay medium. Incubate cells at 37°C with 5% CO₂ overnight.
2. Dilute ionomycin in assay medium to 4 μ M. Prepare threefold serial dilution of PMA in assay medium. Add 25 μ l of diluted ionomycin and 25 μ l of diluted PMA to stimulated wells (Final concentration of ionomycin 1 μ M).
3. Add 50 μ l of assay medium with same concentration of DMSO to the unstimulated control wells (for measuring uninduced level of NF- κ B reporter activity).
4. 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 hours.
6. Add 100 μ l of ONE-Step™ Luciferase Assay reagent per well. Incubate at room temperature for ~15 to 30 minutes and measure luminescence using a luminometer.

Figure 2. PMA dose response in NF- κ B reporter (Luc)-Jurkat cells. Cells were treated with PMA plus ionomycin for ~ 3 hours. The results were shown as fold induction of luciferase

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

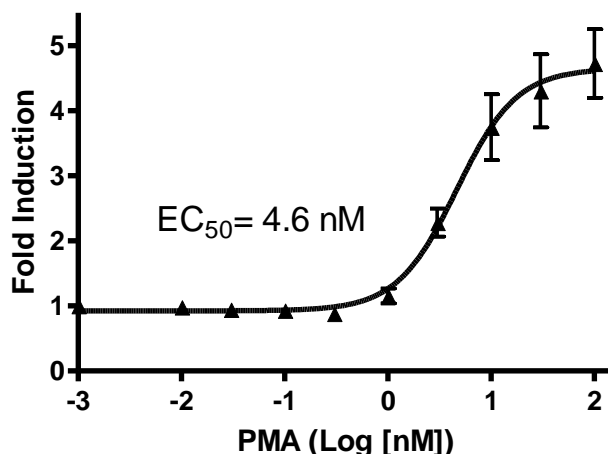
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

reporter expression. Fold induction was determined by comparing values against the mean value for control cells with DMSO treatment.

The EC₅₀ of PMA in the presence of ionomycin in this cell line is 4.6 nM.



Related Products

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
NF-κB reporter (Luc) - HEK293 Cell line	60650	2 vials
NF-κB Reporter (Luc) - A549 Cell Line	60625	2 vials
NF-κB Reporter (Luc) - HCT116 Cell Line	60623	2 vials
NF-κB Reporter (Luc) - CHO-K1 Cell Line	60622	2 vials
CD40/NF-κB Reporter (Luc) - HEK293 Cell Line	60626	2 vials
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
Thaw Medium 2	60184	100 ml
NF-κB Reporter Kit (NF-κB Signaling Pathway)	60614	500 rxns.

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.

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

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