



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

GAS (IFN γ /JAK/STAT1 pathway) Reporter (Luc) – HeLa Cell Line **Catalog #: 79041**

Product description

The GAS reporter (Luc)-HeLa cell line is designed to monitor the activity of interferon gamma-induced signal transduction pathways in cultured cells by measuring activated STAT1 homodimers. It contains a firefly luciferase gene driven by three copies of the interferon gamma-activated sites (GAS) located upstream of the minimal TATA promoter. IFN γ first binds to a heterodimeric receptor consisting of two chains, IFNGR1 and IFNGR2, causing its dimerization and the activation of specific Janus family kinases (JAK1 and JAK2). Two STAT1 molecules associate with this ligand-activated receptor complex and are activated by phosphorylation to form active homodimer. The active STAT1 homodimers translocate to the nucleus where they bind interferon gamma-activated sites (GAS) in the promoter of IFN γ inducible genes, including luciferase reporter gene.

Application

- Monitor interferon gamma-induced signal transduction pathways.
- Screen for activators or inhibitors of JAK/STAT1 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, #MP0025) to confirm the absence of *Mycoplasma* species.

General Culture Conditions

Thaw Medium 1 (BPS Bioscience #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS, 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone #SV30010.01)

Complete Growth Medium: Thaw Medium 1 (BPS Bioscience #60187) and 800 μ g/ml of Geneticin (Life Technologies #11811031).

Cells should be grown at 37°C with 5% CO₂ using complete growth medium.

GAS reporter (Luc)-HeLa 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

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 thaw the cells, it is recommended to quickly thaw the frozen cells directly from liquid nitrogen into a 37°C water-bath, then transfer the entire contents of the vial to a tube containing 10 ml of Thaw Medium 1 (**no Geneticin**). Spin down the cells, remove supernatant and resuspend cells in pre-warmed Thaw Medium 1 (**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 1 (**no Geneticin**). At first passage, switch to complete growth medium (**contains Geneticin**).

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add complete growth medium 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 0.05% Trypsin/EDTA. Add complete growth medium and transfer to a tube, spin down cells, and resuspend in freezing medium (10% DMSO + 90% FBS). Place at -80°C overnight and transfer to 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

- IFN γ (PeproTech #300-02)
- IFN α (PBL Assay Science #11100-1)
- Assay medium: Thaw Medium 1 (BPS Bioscience, #60187) or
- MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS, 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone #SV30010.01)
- 96-well tissue culture treated white clear-bottom assay plate (Corning #3610)
- One-Step luciferase assay system (BPS Bioscience #60690)
- Luminometer

A. IFN γ dose response

1. Harvest GAS reporter (Luc)-HeLa cells from culture in growth medium and seed cells at a density of ~20,000 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 IFN γ in assay medium. Add 50 μ l of diluted IFN γ to IFN γ -stimulated wells.

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

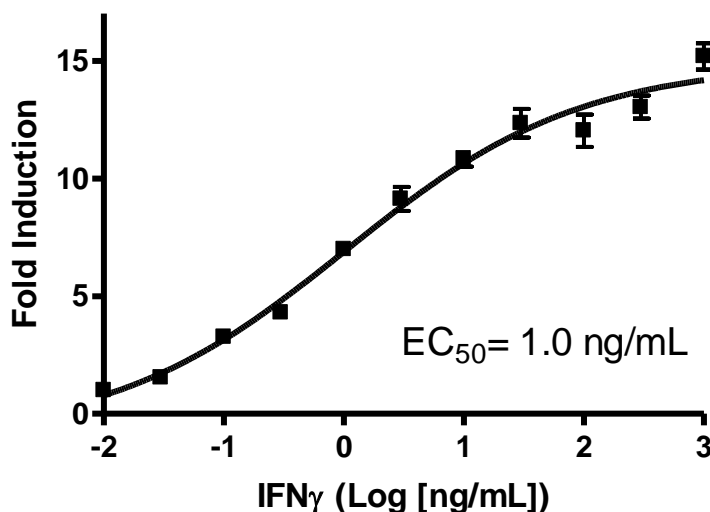
Add 50 μ l of assay medium to the unstimulated control wells (for measuring uninduced level of GAS reporter activity).

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 ~18-24 hours.
6. Prepare ONE-Step™ Luciferase Assay reagent as directed. 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.

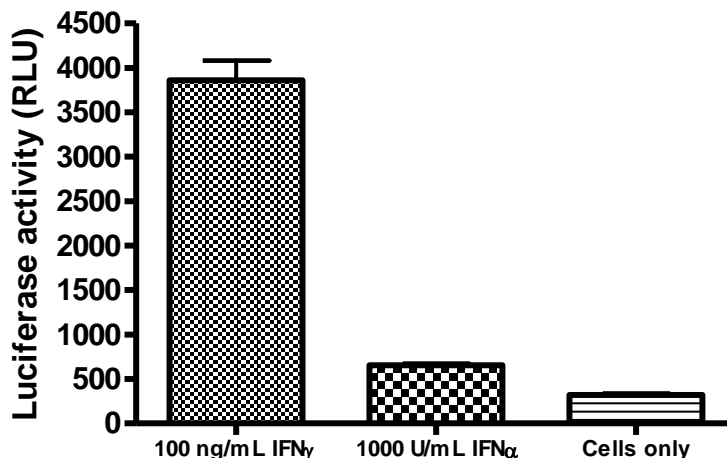
Figure 1. IFN γ dose response in GAS reporter (Luc)-HeLa cells. Cells were treated with IFN γ for ~ 18 hours. The results were shown as fold induction of luciferase reporter expression. Fold induction was determined by comparing values against the mean value for control cells without IFN γ treatment.

The EC₅₀ of IFN γ in this cell line is ~1.0 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

Figure 2. GAS reporter activity in response to IFN α . Cells were seeded at 20,000 cells/well on a white opaque 96-well plate overnight in assay medium before treatment with various human cytokines (IFN γ , 100 ng/ml; IFN α , 1000 ng/ml) and incubated for 18 hours, followed by the addition of luciferin according to manufacturer's protocol (ONE-Step™ Luciferase assay system, BPS Bioscience, #60690-2).



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
Human Interferon-gamma	90162-A	20 μ g
Human Interferon-gamma	90162-B	100 μ g
Mouse Interferon-gamma	90163-A	20 μ g
Mouse Interferon-gamma	90163-B	100 μ g
Human Interferon-alpha 2a	90158-A	20 μ g
Human Interferon-alpha 2b	90159-A	20 μ g
STAT3, GST-tag	75003	20 μ g
ISRE Reporter – HEK293 Recombinant Cell Line	60510	2 vials
ISRE Reporter Kit (JAK/STAT Signaling Pathway)	60613	500 rxns.
Jak1, GST-tag	40449	10 μ g
Jak2 (JH1 domain), His-tag	40450	10 μ g
Jak2 (JH1, JH2 domain), His-tag	40451	10 μ g

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

References

1. Decker T, Kovarik P, Meinke A. (1997) GAS elements: a few nucleotides with a major impact on cytokine-induced gene expression. *J Interferon Cytokine Res.* **17(3)**: 121-34.
2. Darnell J, Kerr IM, Stark GR. (1994) Jak-STAT pathways and transcriptional activation in response to IFNs and other extracellular signaling proteins. *Science* **264**: 1415-1421.

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