



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

PAI-1 Reporter (Luc) – Mv1 Lu Cell line

Catalog #: 60544

Description

PAI-1 Reporter (Luc)-Mv1 Lu cell line is designed for monitoring transforming growth factor β (TGF- β)-induced plasminogen activator inhibitor-1 (PAI-1) expression. Transforming growth factor- β (TGF- β) is a potent regulator of cellular differentiation, proliferation, migration, and protein expression.

PAI-1 Reporter (Luc) –Mv1 Lu cell line contains a firefly luciferase gene under the control of PAI-1 responsive elements stably integrated into Mv1 Lu (NBL-7) cells, showing TGF- β pathway response. This cell line is validated for the TGF- β response to the induction of PAI-1 gene expression through luciferase activity.

Application

- Monitor TGF- β signaling pathway activity

Format

Each vial contains 2×10^6 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 Cat. #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Invitrogen #26140-079), 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 Cat. #60187) and 700 μ g/ml of Geneticin (Invitrogen, #11811031).

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

It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 1 (**no Geneticin**), and spin down the cells. Resuspend cells in pre-warmed Thaw Medium 1 (**no Geneticin**) transfer the resuspended cells

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

To place your order, please contact us by Phone **1.858.829.3082** 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 a T25 flask, and culture in a 37°C CO₂ incubator. At first passage, switch to growth medium containing Geneticin. Cells should be split before they reach complete confluence.

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.25% Trypsin/EDTA, add complete growth medium and transfer to a tube. Spin down cells, resuspend them, and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:10 to 1:20 weekly

Functional Validation and Assay Performance

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

Materials Required but Not Supplied

- Human TGF-β1 (R+D Systems #240-B-002,)
- Thaw Medium 1 (BPS Bioscience, #60187)
- Growth medium: Thaw Medium 1 + G418 700 µg/ml.
- Assay medium: DMEM medium (Hyclone, #SH30243.01) + 1% BSA + 1% Pen/Strep
- 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. Dose response of PAI-1 Reporter (Luc)-Mv1 Lu cells to TGF-β1

1. Harvest PAI-1 Reporter (Luc)-Mv1 Lu cells from culture in Growth medium and seed cells at a density of 20000~30,000 cells per well into white clear-bottom 96-well microplate in 100 µl of growth medium without selection (G418) at 37°C in a CO₂ incubator overnight. Leave a couple of wells empty for the cell free control.

2. Prepare TGF-β1 solution in assay medium (10-point serial dilutions from 100 ng/ml).

3. Remove the growth medium from cells and add 60 µl of diluted TGF-β1 to each test well. Add 60 µl of assay medium to the unstimulated control wells and 60 µl of assay medium to the cell-free control wells (for determining background luminescence). *Set up each treatment in at least triplicate.*

4. Incubate cell plates at 37°C in a CO₂ incubator for ~ 24 hours.

5. Perform luciferase assay using ONE-STEP™ luciferase assay system: Add 60 µl of One-STEP Luciferase reagent per well and rock gently at room temperature for 30 minutes. Measure luminescence using a luminometer.

6. Data Analysis: Subtract average background luminescence (cell-free control wells) from luminescence reading of all 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**

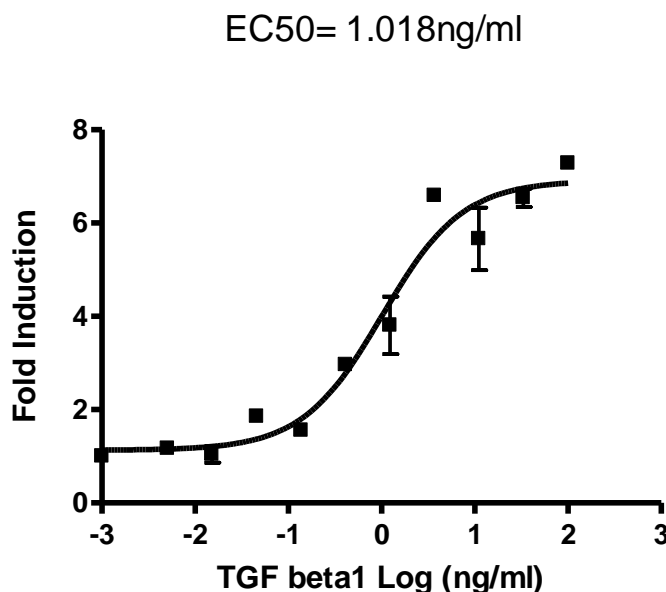
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Fold induction of PAI-1 luciferase reporter expression = background-subtracted luminescence of TGF- β 1-stimulated well / average background-subtracted luminescence of unstimulated control wells

Figure 1. Dose response of PAI-1 Reporter (Luc)-Mv1 Lu cells to human TGF- β 1

The results are shown as fold induction of PAI-1 luciferase reporter expression. The EC50 of TGF- β 1 is ~ 1 ng/ml.



Sequence

PAI-1 promoter sequence (874bp-EcoR1/Hind III)

```
AAGCTTTTACCATGGTAACCCCTGGTCCCGTTCAGCCACCACCACCCACCCAGCACACCT
CCAACCTCAGCCAGACAAGGTTGTTGACACAAGAGAGCCCTCAGGGGCACAGAGAGAGTC
TGGACACGTGGGGAGTCAGCCGTGTATCATCGGAGGCGGCCGGGCACATGGCAGGGATG
AGGAAAGACCAAGAGTCCTCTGTTGGGCCAAGTCCTAGACAGACAAAACCTAGACAAT
CACGTGGCTGGCTGCATGCCCTGTGGCTGTTGGGCTGGGCCAGGAGGAGGGAGGGGC
GCTCTTTCCTGGAGGTGGTCCAGAGCACCGGGTGGACAGCCCTGGGGGAAAACCTCCAC
GTTTTGATGGAGGTTATCTTTGATAACTCCACAGTGACCTGGTTCGCCAAAGGAAAAGCAG
GCAACGTGAGCTGTTTTTTTTTCTCCAAGCTGAACACTAGGGGTCTAGGCTTTTTGGGT
CACCCGGCATGGCAGACAGTCAACCTGGCAGGACATCCGGGAGAGACAGACACAGGCAG
AGGGCAGAAAGGTCAAGGGAGGTTCTCAGGCCAAGGCTATTGGGGTTTGCTCAATTGTTT
CTGAATGCTCTTACACACGTACACACACAGAGCAGCACACACACACACACATGCCTC
AGCAAGTCCCAGAGAGGGAGGTGTCGAGGGGGACCCGCTGGCTGTTTCAAGCGACTCCC
AGAGCCAGTGAGTGGGTGGGGCTGGAACATGAGTTCATCTATTTCTGCCACATCTGGT
ATAAAAGGAGGCAGTGGCCACAGAGGAGCACAGCTGTGTTTGGCTGCAGGGCCAAGAG
CGCTGTCAAGAAGACCCACACGCCCCCTCCAGCAGCTGAATTC
```

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

Abe, M. *et al.* (1994) An assay for transforming growth factor- β using cells transfected with a plasminogen activator inhibitor-1 promoter-luciferase construct. *Analytical Biochemistry* **216**:276-584.

Khan, S. *et al.* (2012) Quantification of active and total transforming growth factor- β levels in serum and solid organ tissues by bioassay. *BMC Research Notes* **5**: 636

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
TGF β 1	90900-1	1 μ g
TGF β 1	90900-2	5 μ g
TGF β 1	90900-10	10 μ g
TGF β 1	90900-3	1 mg
TGF β 1, Latent	90901-1	5 μ g
TGF β 1, Latent	90901-2	25 μ g
TGF β 1, Latent	90901-3	1 mg
TGF β R2, GST-tag	40707	50 μ g
TGF/SMAD Signaling Pathway	60653	2 vials
SBE Reporter – HEK293 Cell Line		

Notes

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