



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

Data Sheet

JNK Signaling Pathway **AP1 Reporter - HEK293 Recombinant Cell line** **Catalog #: 60405**

Description

The stress-activated protein kinase/c-jun N-terminal kinase (SAPK/JNK) family of proteins includes mitogen-activated protein kinases (MAPKs) that are activated by stress, inflammatory cytokines, mitogens, oncogenes, and inducers of cell differentiation and morphogenesis. Upon activation of the SAPK/JNK pathway, MAP Kinase Kinases phosphorylate and activate JNKs. The activated JNKs translocate to the nucleus where they phosphorylate and activate transcription factors such as c-Jun. cJun then binds to the activator protein-1 (AP1) response element and induces AP1 transcription.

The AP1 Reporter – HEK293 cell line contains a firefly luciferase gene under the control of AP1-responsive elements that are stably integrated into HEK293 cells. This cell line is validated for its response to stimulation by Phorbol 12-Myristate 13-Acetate (PMA) and to treatment with inhibitors of the JNK signaling pathway.

Application

- Monitor the JNK signaling pathway activity and AP1-mediated activity.
- Screen for activators or inhibitors of the JNK signaling pathway.

Format

Each vial contains ~1.5 X 10⁶ cells in 1 ml of 10% DMSO.

Storage

Immediately upon receipt, store in liquid nitrogen.

General Culture Conditions

Thaw Medium 1 (BPS Cat. #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Life technologies #26140-079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na-pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone SV30010.01)

Growth Medium 1B (BPS Cat. #79531) Thaw Medium 1 (BPS Cat. #60187) and 400 µg/ml of Geneticin (Life Technologies #11811031).

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 West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 1B (BPS Cat. #79531) (Thaw Medium 1 and Geneticin). It may be necessary to adjust the percentage of CO₂ in the incubator depending on the NaHCO₃ level in the basal medium.

To thaw the cells, 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**). Spin down cells, resuspend cells in pre-warmed Thaw Medium 1 (**no Geneticin**), transfer resuspended cells to a T25 flask and culture in a CO₂ incubator at 37°C. At first passage, switch to Growth Medium 1B (**contains Thaw Medium 1 and 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 using 0.05% Trypsin/EDTA, add Growth Medium 1B transfer to a tube, spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ration: 1:10 to 1:20 weekly.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with Trypsin/EDTA. Add Growth Medium 1B 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.

Functional Validation and Assay Performance

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

Materials Required but Not Supplied for Cell Culture

- Thaw Medium 1 (BPS Cat. #60187)
- Growth Medium 1B (BPS Cat. #79531)

Materials Required but Not Supplied for Cellular Assay

- Phorbol 12-Myristate 13-Acetate (PMA) (LC Laboratories # P-1680). Prepare stock solution in DMSO.
- JNK inhibitor V (AS601245) (Santa Cruz biotechnology # sc-202672): inhibitor of JNK1,2,3. Prepare stock solution in DMSO.
- Assay Medium 1B (BPS Bioscience #79617): Opti-MEM I (Life technologies # 31985-062), 0.5% FBS, 1% non-essential amino acids, 1mM Na pyruvate, and 1% Pen/Strep
- 96-well tissue culture plate or 96-well tissue culture-treated white clear-bottom assay plate
- ONE-Step™ Luciferase Assay Reagent (BPS Cat. #60690)
- Luminometer

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 West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Mycoplasma testing

The cell line has been screened using the PCR-based VenorGeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of Mycoplasma species.

A. Response of AP1 Reporter – HEK293 cells to PMA

1. Harvest AP1 Reporter – HEK293 cells from culture in Growth Medium 1B and seed cells at a density of ~ 35,000 cells per well into a white clear-bottom 96-well microplate in 100 µl of Thaw Medium 1.
2. Incubate cells at 37°C in a CO₂ incubator for overnight.
3. The next day, carefully remove the medium from the wells. Make 3-fold serial dilutions of PMS in Assay Medium 1B. The final concentration of DMSO in assay medium in all dilutions should be 0.1%.

Assay Medium 1B (BPS Bioscience #79617): Opti-MEM I (Life technologies # 31985-062), 0.5% FBS, 1% non-essential amino acids, 1mM Na pyruvate, and 1% Pen/Strep

4. Add 100 µl of serial dilutions of PMA in Assay Medium 1B to stimulated wells. Add 100 µl of Assay Medium 1B with 0.1% DMSO to unstimulated control wells. Add 100 µl of Assay Medium 1B with 0.1% DMSO to cell-free control wells (for determining background luminescence). Set up each treatment in at least triplicate.
5. Incubate the plate at 37°C in a CO₂ incubator for ~ 6 hours.
6. Perform luciferase assay using the ONE-Step™ Luciferase Assay System following the protocol provided: Add 100 µl of ONE-Step™ Luciferase reagent per well and rock at room temperature for ~15 minutes. Measure luminescence using a luminometer.
If using luciferase reagents from other vendors, follow the manufacturer's assay protocol.
7. Data Analysis: Subtract average background luminescence (cell-free control wells) from luminescence reading of all wells.
The fold induction of AP1 luciferase reporter expression = background-subtracted luminescence of stimulated well / average background-subtracted luminescence of unstimulated control wells

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

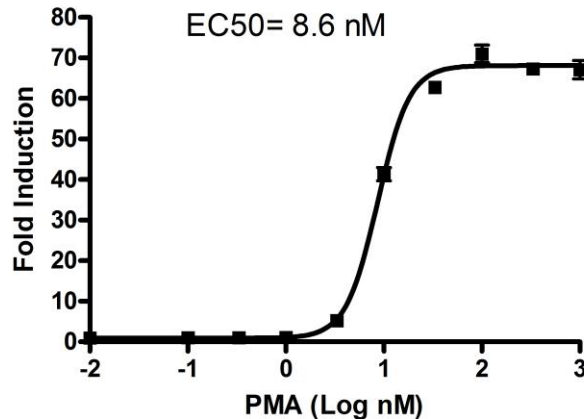
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

Figure 1. Dose response of AP1 Reporter – HEK293 cells to PMA.

The results are shown as fold induction of AP1 luciferase reporter expression.

**B. Inhibition of PMA-induced reporter activity by an inhibitor of the JNK signaling pathway in AP1 Reporter – HEK293 cells**

1. Harvest AP1 Reporter – HEK293 cells from culture in Growth Medium 1B and seed cells at a density of 35,000 cells per well into a white clear-bottom 96-well microplate in 100 μ l of Thaw Medium 1.
2. Incubate cells at 37°C in a CO₂ incubator for overnight.
3. The next day, dilute the inhibitor (JNK inhibitor V) stock in Assay Medium 1B. Carefully remove the medium from wells and add 90 μ l of diluted inhibitor in Assay Medium 1B to the wells. The final concentration of DMSO in Assay Medium 1B can be up to 0.5%.

Assay Medium 1B (BPS Bioscience #79617): Opti-MEM I (Life technologies # 31985-062), 0.5% FBS, 1% non-essential amino acids, 1mM Na pyruvate, and 1% Pen/Strep

Add 90 μ l of Assay Medium 1B with same concentration of DMSO without inhibitor to inhibitor control wells.

Add 90 μ l of Assay Medium 1B with DMSO to cell-free control wells (for determining background luminescence).

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 West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

4. Incubate the plate at 37°C in a CO₂ incubator for 1 hour.
5. Add 10 µl of diluted PMA in Assay Medium 1B to stimulated wells (final [PMA] = 10 nM). Final DMSO concentration should be 0.1%.
Add 10 µl of Assay Medium 1B with 0.1% DMSO to the unstimulated control wells (cells without inhibitor and PMA treatment for determining the basal activity).
Add 10 µl of Assay Medium 1B with 0.1% DMSO to cell-free control wells.
Set up each treatment in at least triplicate.

Treatment Reference Guide

	Stimulated Wells		Unstimulated Control Wells	Cell-free Control Wells
	With inhibitor	Without inhibitor (control well)		
Step 3	90 µl diluted inhibitor in Assay Medium 1B	90 µl Assay Medium 1B with DMSO only	90 µl Assay Medium 1B with DMSO only	90 µl Assay Medium 1B with DMSO only
Step 5	10 µl PMA in Assay Medium 1B (final [PMA] = 10 nM)	10 µl PMA in Assay Medium 1B (final [PMA] = 10 nM)	10 µl Assay Medium 1B with 0.1% DMSO	10 µl Assay Medium 1B with 0.1% DMSO

6. Incubate the plate at 37°C in a CO₂ incubator for ~6 hours.
7. Perform luciferase assay using the ONE-Step™ Luciferase Assay System following the protocol provided: Add 100 µl of ONE-Step™ Luciferase reagent per well and rock at room temperature for ~15 minutes. Measure luminescence using a luminometer.
If using other luciferase reagents from other vendors, follow the manufacturer's assay protocol.
8. Data Analysis: Obtain background-subtracted luminescence by subtracting 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.

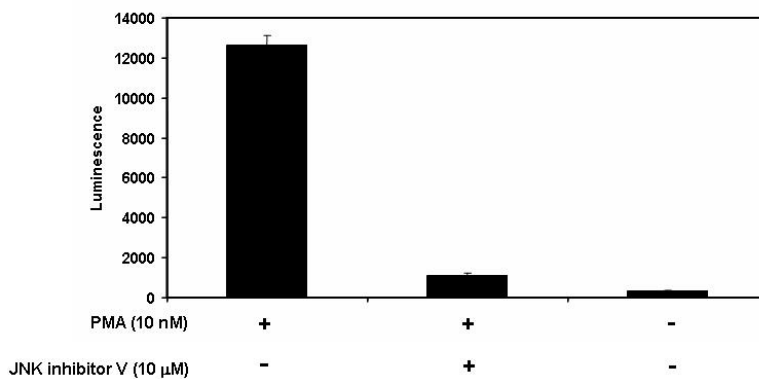
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

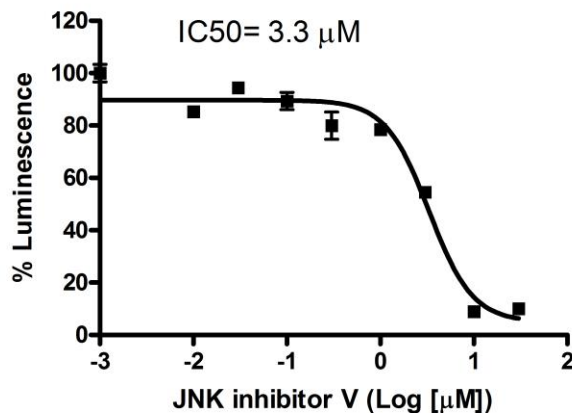
Figure 2. Inhibition of PMA-induced reporter activity by a JNK pathway inhibitor in AP1 Reporter – HEK293 cells

2a. JNK inhibitor V blocked PMA-induced AP1 reporter activity.



2b. JNK inhibitor V inhibition dose response curve

The results are shown as percentage of luminescence. The background-subtracted luminescence of cells stimulated with PMA in the absence of JNK inhibitor V was set at 100%.



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 West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

References:

1. Zhou H. *et. al.* (2005) Frequency and distribution of AP-1 sites in the human genome. *DNA Research*. **11**: 139-150.
2. Gaillard P. *et.al.* (2005) Design and synthesis of the first generation of novel potent, selective, and in vivo active (benzothiazol-2-yl)acetonitrile inhibitors of the c-Jun N-terminal kinase. *J Med Chem*. **48(14)**:4596-4607.

Related Products

<u>Product Name:</u>	<u>Catalog #</u>	<u>Size</u>
Thaw Medium 1	60187	100ml
Growth Medium 1B	79531	500 ml
AP1 Reporter Kit (JNK Signaling Pathway)	60612	500 reactions
SRE Reporter - HEK293 Cell line (ERK Pathway)	60406	2 vials
SRE Reporter Kit (MAPK/ERK Signaling Pathway)	60511	500 reactions
MAPK10 (JNK3), human	40092	10 µg
JNK1-β1(K55M), human	40871	100 µg
MAP3K14 (NIK), human	40090	10 µg
MAPKAPK2 (MK2), human	40088	100 µg
JNK1, mouse	40071	10 µg
JNK2, human	40113	10 µg
JNK3, human	40114	10 µg
ERK1, human	40055	10 µg
ERK2, human	40299	10 µg
ERK2, inactive, human	40056	10 µg
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml

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