



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

AP1 Reporter Kit (JNK Signaling Pathway) **Catalog #: 60612**

Background

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. The activated c-Jun forms homodimers or heterodimers with fos family proteins which bind to the activator protein-1 (AP1) response element and induce target gene transcription.

Description

The *AP1 Reporter Kit* is designed for monitoring the activity of the JNK signaling pathway and the transcriptional activity of AP1 in cultured cells. The kit contains a transfection-ready AP1 luciferase reporter vector. This reporter contains the firefly luciferase gene under the control of multimerized AP1 responsive elements located upstream of a minimal promoter. The AP1 reporter is premixed with a constitutively-expressing *Renilla* luciferase vector that serves as an internal control for transfection efficiency.

The kit also includes a non-inducible firefly luciferase vector premixed with constitutively-expressing *Renilla* luciferase vector as a negative control. The non-inducible luciferase vector contains the firefly luciferase gene under the control of a minimal promoter, without any additional response elements. The negative control is critical for determining pathway-specific effects and the background luciferase activity.

For assays in more difficult to transfect cell types, we suggest using our lentiviral products. The corresponding lentiviral product for this kit is BPS Bioscience #79823.

Applications

- Monitor JNK signaling pathway activity and AP1-mediated activity.
- Screen for activators or inhibitors of the JNK signaling pathway.
- Study effects of RNAi or gene overexpression on the activity of the JNK pathway.

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Components

Component	Specification	Amount	Storage
Reporter (Component A)	AP1 luciferase reporter vector + constitutively expressing <i>Renilla</i> luciferase vector	500 µl (60 ng DNA/ µl)	-20°C
Negative Control Reporter (Component B)	Non-inducible luciferase vector + constitutively expressing <i>Renilla</i> luciferase vector	500 µl (60 ng DNA/ µl)	-20°C

Note: These vectors are designed for transient transfection. They are NOT SUITABLE for transformation and amplification in bacteria.

Materials Required but Not Supplied

- Mammalian cell line and appropriate cell culture medium
- 96-well tissue culture plate or 96-well tissue culture-treated white clear-bottom assay plate
- Transfection reagent for mammalian cell line [We use Lipofectamine™ 2000 (Life Technologies #11668027). However, other transfection reagents work equally well.]
- Opti-MEM I Reduced Serum Medium (Life Technologies #31985-062)
- TWO-step luciferase assay:
 TWO-Step Luciferase (Firefly & Renilla) Assay System (BPS Bioscience #60683):
 This system assays cells directly in growth medium. It can be used with any luminometer. Automated injectors are not required.
- Luminometer

Generalized Transfection and Assay Protocols

The following procedure is designed to transfect the reporter into HEK293 or HeLa cells using Lipofectamine 2000 in a 96-well format. To transfect cells in different tissue culture formats, adjust the amounts of reagents and cell number in proportion to the relative surface area. If using a transfection reagent other than Lipofectamine 2000, follow the manufacturer's transfection protocol. Transfection conditions should be optimized according to the cell type and study requirements.

All amounts and volumes in the following setup are given on a per-well basis.

1. One day before transfection, seed cells at a density of ~ 30,000 cells per well in 100 µl of growth medium so that cells will be 90% confluent at the time of transfection.
2. The next day, for each well, prepare complexes as follows:
 - a. Dilute DNA mixtures in 15 µl of Opti-MEM I medium (antibiotic-free). Mix gently.

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Depending upon the experimental design, the DNA mixtures may be any of following combinations:

- **1 µl of Reporter** (component A); in this experiment, the control transfection is **1 µl of Negative Control Reporter** (component B).
- **1 µl of Reporter** (component A) + experimental vector expressing gene of interest; in this experiment, the control transfections are: **1 µl of Reporter** (component A) + negative control expression vector, **1 µl of Negative Control Reporter** (component B) + experimental vector expressing gene of interest, and **1 µl of Negative Control Reporter** (component B) + negative control expression vector.
- **1 µl of Reporter** (component A) + specific siRNA; in this experiment, the control transfections are: **1 µl of Reporter** (component A) + negative control siRNA, **1 µl of Negative Control Reporter** (component B) + specific siRNA, and **1 µl of Negative Control Reporter** (component B) + negative control siRNA.

Note: we recommend setting up each condition in at least triplicate and preparing transfection cocktail for multiple wells to minimize pipetting errors.

- b. Mix Lipofectamine 2000 gently before use, then dilute 0.35 µl of Lipofectamine 2000 in 15 µl of Opti-MEM I medium (antibiotic-free). Incubate for 5 minutes at room temperature.

Note: Prepare this dilution cocktail in volumes sufficient for the whole experiment.

- c. After the 5 minute incubation, combine the diluted DNA with diluted Lipofectamine 2000. Mix gently and incubate for 25 minutes at room temperature.

3. Add the 30 µl of the complexes to each well containing cells and medium. Mix gently by tapping the plate.
4. Incubate cells at 37°C in a CO₂ incubator for overnight.
5. The next day, change medium to assay medium (Opti-MEM I, 0.5% FBS, 1% non-essential amino acids, 1 mM Na-pyruvate, 1% Pen/Strep) containing an activator of AP1 such as PMA. Incubate cells at 37°C in a CO₂ incubator for ~ 6 to 24 hours. After treatment, perform the TWO-Step Luciferase (Firefly & Renilla) Assay System following the protocol on the BPS data sheet (BPS Bioscience #60683).

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Sample protocol to determine the effect of PMA on AP1 reporter activity in HEK293 or HeLa cells

1. One day before transfection, seed HEK293 or HeLa cells at a density of 30,000 cells per well into white clear-bottom 96-well plate in 100 μ l of growth medium (MEM/EBSS (Hyclone #SH30024.01), 10% FBS, 1% non-essential amino acids, 1 mM Na-pyruvate, 1% Pen/Strep). Incubate cells overnight at 37°C in a CO₂ incubator.
2. The next day, transfect 1 μ l of AP1 reporter (component A) into cells following the procedure in **Generalized Transfection and Assay Protocols**.
3. Incubate cells at 37°C in a CO₂ incubator overnight.
4. The next day after transfection, prepare 1 mM stock solution of PMA in DMSO. Dilute the PMA stock in assay medium. Change cell medium to 50 μ l of diluted PMA in assay medium to induce the AP1 reporter. For unstimulated control wells, treat cells with 50 μ l of assay medium (without PMA). Add 50 μ l of assay medium (without PMA) to cell-free control wells to determine the background luminescence. Set up each treatment in at least triplicate.
5. Incubate cells at 37°C in a CO₂ incubator for ~6 hours.
6. Perform TWO-step luciferase assay using TWO-Step Luciferase (Firefly & Renilla) Assay System (BPS Bioscience #60683): Dilute 100x Firefly Luciferase Reagent Substrate (Component B) into Firefly Luciferase Reagent Buffer (Component A). Add 50 μ l of Firefly Luciferase reagent per well and rock at room temperature for ~15 minutes, then measure firefly luminescence using a luminometer. Dilute 100x Renilla Luciferase Reagent Substrate (Component D) into Renilla Luciferase Reagent Buffer (Component C). Add 50 μ l of Renilla Luciferase reagent per well, rock at room temperature for ~1 minute and measure Renilla luminescence.
7. To obtain the normalized luciferase activity for the AP1 reporter, subtract the background luminescence, then calculate the ratio of firefly luminescence from AP1 reporter to *Renilla* luminescence from the control *Renilla* luciferase vector.

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

Figure 1. PMA induced the expression of AP1 reporter. The results are shown as fold induction of normalized AP1 reporter activity. Fold induction is determined by comparing values against the mean value for unstimulated control cells.

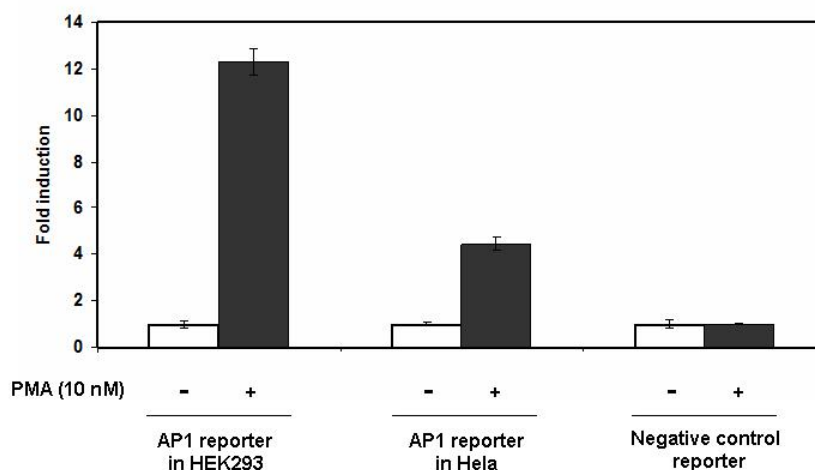
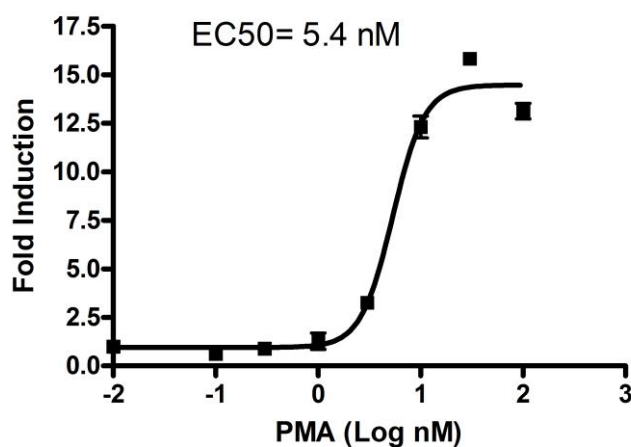


Figure 2. Dose response of AP1 reporter activity to PMA in HEK293 cells. The results are shown as fold induction of normalized AP1 reporter activity. Fold induction is determined by comparing values against the mean value for unstimulated control cells. The EC₅₀ of PMA is ~5.4 nM



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: support@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

Sample protocol to determine the effect of JNK pathway inhibitor on AP1 reporter activity

1. One day before transfection, seed cells at a density of 30,000 cells per well into white clear-bottom 96-well plate in 100 μ l of growth medium. Incubate cells overnight at 37°C in a CO₂ incubator.
2. The next day, transfect 1 μ l of AP1 reporter (component A) into cells following the procedure in **Generalized Transfection and Assay Protocols**.
3. Incubate cells at 37°C in a CO₂ incubator for overnight.
4. The next day after transfection, prepare 10 mM stock solution of JNK inhibitor V in DMSO. Dilute the inhibitor in assay medium to a final concentration of 10 μ M. Carefully remove the medium from wells and add 45 μ l of diluted inhibitor in assay medium to the wells. Add 45 μ l of assay medium (without inhibitor) to inhibitor control wells or unstimulated control wells. Add 45 μ l of assay medium (without inhibitor) to cell-free control wells (for determining background luminescence). Set up each treatment in at least triplicate.
5. Incubate cells at 37°C in a CO₂ incubator for 1 hour.
6. Add 5 μ l of diluted PMA in assay medium to wells. The final PMA concentration for the cells is 10 nM. Add 5 μ l of assay medium without PMA to the unstimulated control wells (cells without inhibitor and without PMA treatment for determining the basal activity). Add 5 μ l of assay medium without PMA to cell-free control wells.
7. Incubate cells at 37°C in a CO₂ incubator for 6 hour.
8. Perform TWO-step luciferase assay using TWO-Step Luciferase (Firefly & Renilla) Assay System (BPS Bioscience #60683): Dilute 100x Firefly Luciferase Reagent Substrate (Component B) into Firefly Luciferase Reagent Buffer (Component A). Add 50 μ l of Firefly Luciferase reagent per well and rock at room temperature for ~15 minutes, then measure firefly luminescence using a luminometer. Dilute 100x Renilla Luciferase Reagent Substrate (Component D) into Renilla Luciferase Reagent Buffer (Component C). Add 50 μ l of Renilla Luciferase reagent per well, rock at room temperature for ~1 minute and measure Renilla luminescence.
9. To obtain the normalized luciferase activity for the AP1 reporter, subtract the background luminescence, then calculate the ratio of firefly luminescence from the AP1 reporter to *Renilla* luminescence from the control *Renilla* luciferase vector.

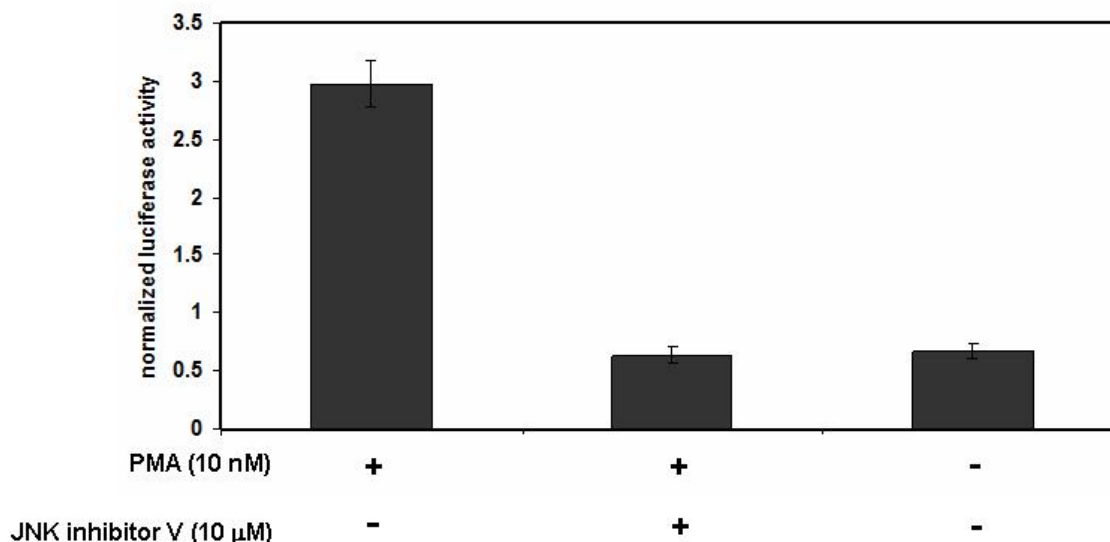
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com

Figure 3. Inhibition of PMA-induced AP1 reporter activity by JNK inhibitor V. The results are shown as normalized AP1 reporter luciferase activity.



References

Zhou H. *et. al.* (2005) Frequency and distribution of AP-1 sites in the human genome. *DNA Research*. **11**: 139-150.

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>
AP1 Reporter – HEK 293 cell line	60405	1 vial
SRE Reporter Kit (MAPK/ERK Signaling Pathway)	60511	500 reactions
TWO-Step Luciferase System (Firefly & Renilla) Assay	60683-1	10 mL
TWO-Step Luciferase System (Firefly & Renilla) Assay	60683-2	100 mL
TWO-Step Luciferase System (Firefly & Renilla) Assay	60683-3	1 L
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

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com