



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

cAMP/PKA Signaling Pathway CRE/CREB Reporter (Luc) – HEK293 Cell Line Catalog #: 60515

Background

The *cAMP/PKA Signaling Pathway CRE/CREB Reporter (Luc) – HEK293 Cell Line* is designed for monitoring the activity of the cAMP/ PKA signaling pathway. The main role of the cAMP response element, or CRE, is mediating the effects of Protein Kinase A (PKA) by way of transcription. It is the main binding site of cAMP response element binding protein (CREB) and is responsible for its activation. CRE is the target of many extracellular and intracellular signaling pathways, including cAMP, calcium, GPCR (G-protein coupled receptors) and neurotrophins. The cAMP/PKA signaling pathway is critical to numerous life processes in living organisms. In the cAMP/PKA signaling pathway, CREB is activated via phosphorylation of PKA and binds to CRE with a general motif of 5'-TGACGTCA-3'. Since CRE is a modulator of the cAMP/PKA signaling pathway, it allows the effects of various inhibitors to be studied.

Description

The *cAMP/PKA Signaling Pathway CRE/CREB Reporter (Luc) – HEK293 Cell Line* contains a firefly luciferase gene under the control of multimerized cAMP response element (CRE) stably integrated into HEK293 cells. Elevation of the intracellular cAMP level activates cAMP response element binding protein (CREB) to bind CRE and induces the expression of luciferase. This cell line is validated for response to stimulation by Forskolin and to the treatment with an inhibitor of the cAMP/PKA signaling pathway.

Application

- Monitor cAMP/PKA signaling pathway activity.
- Screen activators or inhibitors of cAMP/PKA signaling pathway.

Format

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

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.

Storage

Immediately upon receipt, store in liquid nitrogen.

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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

General Culture conditions

Cells should be grown at 37°C with 7% CO₂ using 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), and 100 µg/ml of Hygromycin B (Hyclone #SV30070.01). If culturing cells in medium from other vendors, it may be required to lower the percentage of CO₂ in the incubator depending on the NaHCO₃ level in the basal 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 HEK293 Cell Thaw Medium (BPS Cat #60187), spin down cells (1000 rpm), and resuspend cells in pre-warmed HEK293 Cell Thaw Medium (BPS Cat #60187). Transfer resuspended cells to a single T25 flask and culture at 37°C in a 7% CO₂ incubator. At first passage, switch to HEK293 Cell Thaw Medium (BPS Cat #60187) containing Hygromycin. 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 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.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 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 place in liquid nitrogen the next day. Alternatively, vials may be placed directly in liquid nitrogen.

Note: Just after thawing and at low density, the cells may grow at a slower rate. It is recommended to split the cells with ~ 1:4 ratio at the beginning of culturing. After several passages, the cell growth rate increases and the cells can be split with 1:8 -1:20 ratio weekly.

Functional Validation and Assay Performance

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

Materials Required but Not Supplied

- HEK293 Cell Thaw Medium (BPS Cat #60187)
- Forskolin (LC Laboratories #F-9929), 10 mM in DMSO: activator of cAMP expression
- H-89 (Enzo Life Sciences # BML-E1196-0005) 10 mM in DMSO: inhibitor of cAMP/PKA pathway
- Assay medium: MEM medium (Hyclone #SH30024.01) + 10% FBS + 1% non-essential amino acids + 1 mM Na pyruvate + 1% Pen/Strep
- 96-well tissue culture treated white clear-bottom assay plate (Corning #3610)
- ONE-Step™ Luciferase Assay System (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

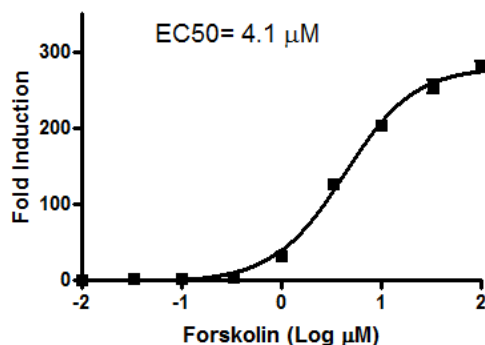


6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

A. Dose response of *cAMP/ PKA Signaling Pathway CRE/CREB Reporter (Luc)* – HEK293 cells to Forskolin

1. Harvest CRE/ CREB reporter (Luc)-HEK293 cells from culture in growth medium and seed cells at a density of ~30,000 cells per well in 45 μ l of assay medium into a white clear-bottom 96-well microplate.
2. Incubate cells at 37°C in a CO₂ incubator overnight (~ 16-18 hours).
3. Add 5 μ l of threefold serial dilution of Forskolin in assay medium to stimulated wells. Add 5 μ l of assay medium to the unstimulated control wells. Add 50 μ l of assay medium to cell-free control wells (for determining background luminescence). Set up each treatment in at least triplicate.
4. Incubate the plate at 37°C in a CO₂ incubator for 5-6 hours.
5. Perform luciferase assay using ONE-Step™ Luciferase Assay System according to the protocol provided: Add 100 μ l of ONE-Step™ Luciferase reagent per well and rock at room temperature for ~10 minutes. Measure luminescence using a luminometer. *If using other luciferase reagents from other vendors, follow the manufacturer's assay protocol.*
6. Data Analysis: Subtract average background luminescence (cell-free control wells) from luminescence reading of all wells.
Fold induction of CRE/CREB luciferase reporter expression = background-subtracted luminescence of Forskolin-stimulated wells / average background-subtracted luminescence of unstimulated control wells

Figure 1. Dose response of CRE/CREB reporter (luc)-HEK293 cells to Forskolin.
The results are shown as fold induction of CRE/CREB 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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

B. Inhibition of Forskolin-induced reporter activity by an inhibitor of cAMP/ PKA signaling pathway in *cAMP/PKA Signaling Pathway CRE/CREB reporter (Luc)-HEK293 cells*

1. Harvest CRE/CREB reporter (Luc)-HEK293 cells from culture in Growth medium and seed cells at a density of ~30,000 cells per well in 40 μ l of assay medium into a white clear-bottom 96-well microplate.
2. Add 5 μ l of inhibitor in assay medium to inhibited wells.
Add 5 μ l of assay medium to the uninhibited control wells.
Set up each treatment in at least triplicate.

Incubate cells at 37°C in a CO₂ incubator overnight (~ 16-18 hours).

3. Add 5 μ l of diluted Forskolin in assay medium to stimulated wells (final Forskolin concentration = 10 μ M).
Add 5 μ l of assay medium to the unstimulated control wells.
Add 50 μ l of assay medium to cell-free control wells (for determining background luminescence).
Set up each treatment in at least triplicate.
4. Incubate the plate at 37°C in a CO₂ incubator for 5-6 hours.
5. Perform luciferase assay using ONE-Step™ Luciferase Assay System according to the protocol provided: Add 100 μ l of ONE-Step™ Luciferase reagent per well and rock at room temperature for ~10 minutes. Measure luminescence using a luminometer.
If using other luciferase reagents from other vendors, follow the manufacturer's assay protocol.
6. 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**

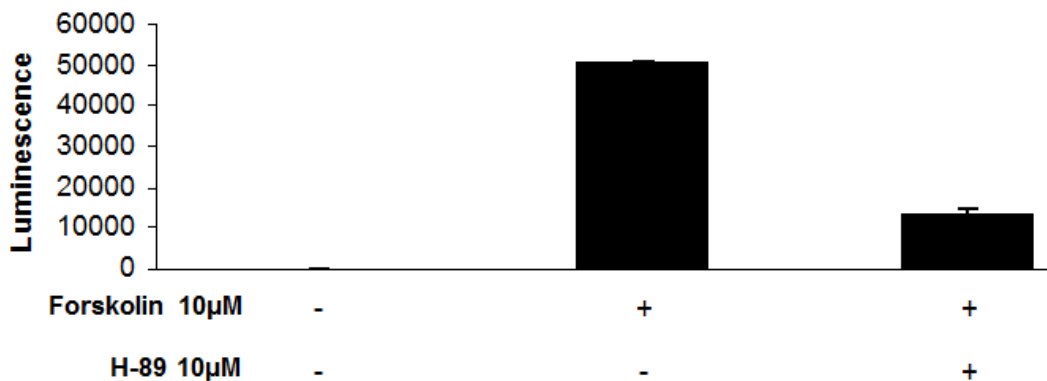
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Figure 2. Inhibition of Forskolin-induced reporter activity by H-89 in CRE/CREB reporter (Luc)-HEK293 cells



Reference(s)

1. Montminy, M.R. *et al.* (1987) Binding of a nuclear protein to the cyclic-AMP response element of the somatostatin gene. *Nature* **328(6126)**: 175-178.
2. Fan Chung, K. (2006) Phosphodiesterase inhibitors in airways disease. *Eur. J. Pharmacol.* **533(1-3)**: 110-117.
3. Malik, R. *et al.* (2008) Cloning, stable expression of human phosphodiesterase 7A and development of an assay for screening of PDE7 selective inhibitors. *Appl. Microbiol. Biotechnol.* **77 (5)**: 1167-1173.

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