



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

PDE10A-HEK293 Recombinant Cell Line

Catalog #: 60410

Lot #: 161121

Description

Recombinant HEK293 cell line expressing human PDE10A2 (phosphodiesterase 10A, accession number NM_006661).

Format

Each vial contains 1×10^6 cells in 1 ml of 10% DMSO.

Storage

Store cells in liquid nitrogen upon arrival. Avoid freeze/thaw cycles.

Introduction

PDE10A plays a key role in signal transduction by regulating the intracellular concentration of cyclic nucleotides. PDE10A hydrolyzes both cAMP and cGMP to nucleoside 5'-monophosphate, but exhibits higher affinity for cAMP, and it utilizes cAMP more efficiently as a substrate (Fujishige K. *et al.*, 1999).

cAMP is an critical second messenger that is required for the proper biological responses of cells to hormones and other extracellular signals. It is required for cellular communication in the hypothalamus/pituitary gland axis and for the feedback control of hormones (Alewijnse A.E. *et al.*, 1997). cAMP is also involved in the activation of protein kinases and ion channel regulation.

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) plus 400 µg/ml of Geneticin (Invitrogen) to ensure maintenance of recombinant PDE10A.

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

Cells should be grown at 37° with 7% CO₂ using complete growth medium. PDE10A-HEK293 cells exhibit a typical cell division time of 24 hours.

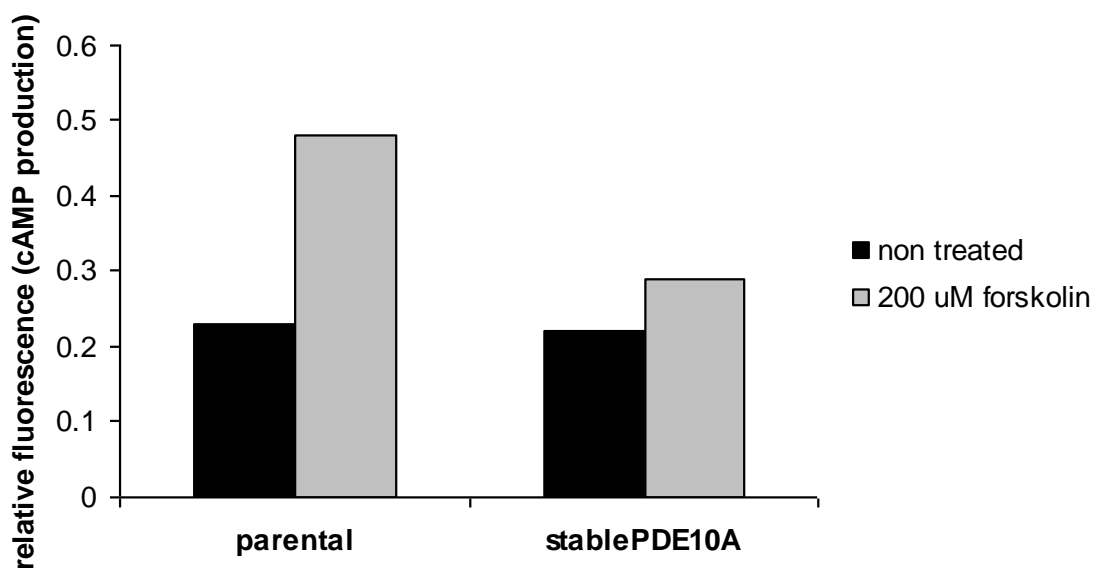
It is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water bath, transfer to a tube containing 10 ml of complete growth medium, spin down cells, resuspend cells and transfer to a T25 flask. Cells should be split prior to reaching complete confluence.

Functional Validation

N-terminal His-tagged human PDE10A2 has been stably expressed in a human embryonic kidney (HEK293) cell line. PDE10A2 expression was confirmed by Western blotting. The function of PDE10A2 was confirmed by cAMP detection assay.

Forskolin is commonly used to raise levels of cyclic AMP (cAMP) in the study and research of cell physiology. Forskolin re-sensitizes cell receptors by activating the enzyme adenylyl cyclase and increasing intracellular levels of cAMP. When HEK293 cells are activated by forskolin, cAMP levels are up-regulated in parental cells whereas cells overexpressing PDE10A show reduction in cAMP accumulation. Inhibition of PDE10A activity restores cAMP levels.

The data provided below demonstrates the effect of PDE10A2 overexpression on cAMP accumulation following forskolin stimulation in HEK293 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

Figure 1. PDE10A2 overexpression reduces cAMP accumulation following forskolin stimulation.

Cells were incubated in PBS buffer for 15 min. and subsequently stimulated with 200 μ M of forskolin for 30 min. cAMP production was measured by detection of fluorescent quenching using a cAMP fluorescent assay kit (Mediomics, St. Louis, MO).

Addition of the pan-PDE inhibitor IBMX to the assay buffer reverses the effect of PDE10A overexpression.

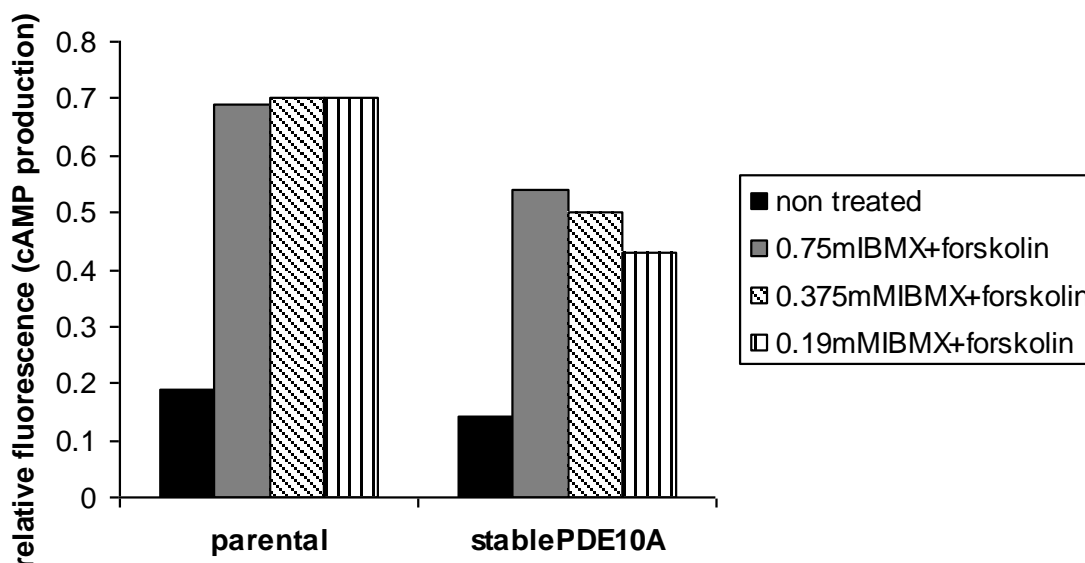


Figure 2. Presence of IBMX, a pan-PDE inhibitor, affects PDE10A2 activity and restores accumulation of cAMP.

Cells were incubated in assay buffer with different concentration of IBMX for 15 min and consequently stimulated with 200 μ M of forskolin for 30 min. cAMP production was measured by detection of fluorescent quenching (Mediomics).

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

Vector and Sequence

Human His-PDE10A2 cloned in pcDNA3.1 (accession number NM_006661)

His-PDE10A sequence:

MHHHHHHHRIEERKSQHLTGLTDEKVKAYLSLHPQVLDEFVSESVSAETVEKWLKRKN
NKSEDESAPKEVSRYQDTNMQGVVYELNSYIEQRLDTGGDNQLLLYELSSIIKIATKAD
GFALYFLGECNNSLCIFTTPPGIKEGKPRIPAGPITQGTTVSAYVAKSRKTLLEDILGDE
RFPRGTGLESGETRIQSVLCLPIVTAIGDLIGILELYRHWGKEAFCLSHQEVATANLAWAS
VAIHQVQVCRGLAKQTELNDFLLDVSKTYFDNIVAIDSLLEHIMIYAKNLVNADRCALFQ
VDHKNKELYSDLFDIGEEKEGKPVFKKTKEIRFSIEKGIAGQVARTGEVLNIPDAYADPR
FNREVDLYTGYTTRNLCMPIVSRGSGVIGVVQMVNKGSGSAFSKTDENNFKMFAVFCAL
ALHCANMYHRIRHSECIYRVTMEKLSYHSICTSEEWQGLMQFTLPVRLCKEIELFHFDI
GPFENMWPGIFVYVMVHRSCGTSCFELEKLCRFIMSVKKNYRRVPYHNWKHAVTVAHC
MYAILQNNHTLFTDLERKGLLIACLDLHRGFSNSYLQKFDHPLAALYSTSTMEQHH
FSQTVSILQLEGHNIFSTLSSSEYEQVLEIRKAIATDLALYFGNRKQLEEMYQTGSLNL
NNQSHRDRVIGLMMTACDLCSVTKLWPVTKLTANDIYAEEFWAEGDEMCKLGIQPIPM
DRDKKDEVPPQGLGFYNAVAIPCYTTLTQILPPTPELLKACRDNLSQWEKVIRGEETAT
WISSPSVAQKAAASED.

References

1. Fujishige K, *et al.* (1999). "Cloning and characterization of a novel human phosphodiesterase that hydrolyzes both cAMP and cGMP (PDE10A)." *J. Biol. Chem.* **274** (26): 18438–45.
2. Alewijnse, AE, *et al.* (1997). "Modulation of forskolin-mediated adenylyl cyclase activation by constitutively active G_s-coupled receptors." *FEBS Letters* **419** (2-3): 171-174.

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

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