



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

Data Sheet

CFTR – HEK293 Recombinant Cell Line

Catalog #: 60506

Background

Cystic fibrosis transmembrane conductance regulator (CFTR) is a protein that in humans is encoded by the CFTR gene. CFTR is an ABC transporter-class ion channel that transports chloride and thiocyanate ions across epithelial cell membranes. Mutations of the CFTR gene affect functioning of the chloride ion channels in these cell membranes, leading to cystic fibrosis [1]. It is characterized by the triad of chronic bronchopulmonary disease (with recurrent respiratory infections), pancreatic insufficiency (which leads to malabsorption and growth retardation) and elevated sweat electrolytes [2].

Description

The CFTR-HEK293 Cell Line expresses full length, human cystic fibrosis transmembrane conductance regulator (CFTR) protein (Genbank # P13569). The expression of CFTR is confirmed by Western blotting.

Sequence

A synthetic codon-optimized DNA sequence encoding human CFTR protein [2] with C-terminal Streptavidin-Binding Peptide (SBP) [3] tag is stably integrated in 293HEK-Trex cells.

Application

- Drug compound screening
- Functional Assays
- Efficient antigen for mouse immunization

Format

Each vial contains 2×10^6 cells in 1 ml of Sigma Freezing Medium (Cat. # C-6164).

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**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)



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

Host cell

293HEK-Trex

Recommended Storage

Immediately upon receipt, store in liquid nitrogen.

Propagation Medium

Growth Medium 4A (BPS Cat. #79535):

Thaw Medium 4 (BPS Cat. #60181) + 10 µg/ml Blasticydin, 0.2 mg/ml Zeocin

Thaw Medium 4, Blasticydin, Zeocin: DMEM/F12 (1:1) (Hyclone # SH30271.01) supplemented with 10% FBS (Life Technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone # SV30010) and 10 µg/ml Blasticydin, 0.2 mg/ml Zeocin.

Induction of target protein expression

Growth Medium 4C (BPS Cat. #79631):

Thaw Medium 4 (BPS Cat. #60181) + 1 µg/ml Doxycycline (Biochemika #44577) and 3 mM Na-butyrate (Acros Organics #263190250):

Thaw Medium 4, Doxycycline, Na-butyrate: DMEM/F12 (1:1) (Hyclone # SH30271.01) supplemented with 10% FBS (Life Technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone # SV30010) 1 µg/ml Doxycycline (Biochemika #44577) and 3 mM Na-butyrate (Acros Organics #263190250)

During 20-24 hrs. before cell harvesting.

Stability

Stable after minimum of seven continuous passages. Upon receipt, amplify the cells in culture and make several frozen aliquots for future use.

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**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)

Figure 1. Western Blot of the CFTR expressing cells. 1. CFTR-GFP-SBP fusion, 2. CFTR-SBP fusion; detection via Streptavidin-AP.

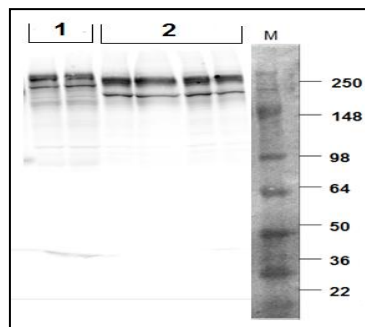
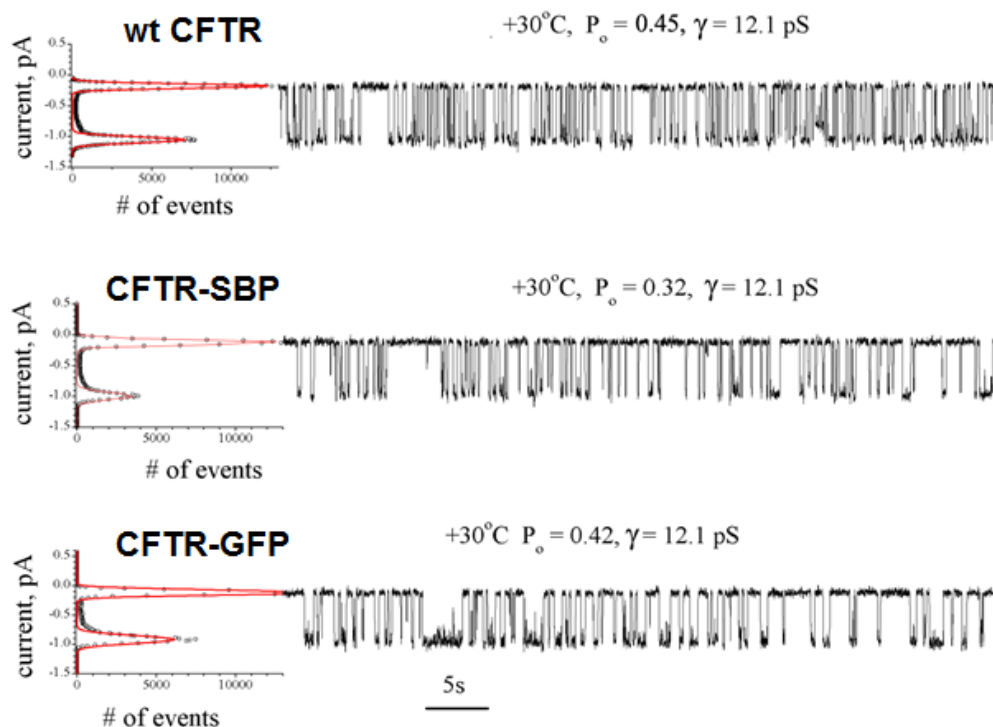


Figure 2. Electrophysiological analysis of CFTR cells. Comparison of wt CFTR vs. CFTR-SBP and CFTR-GFP.



OUR F

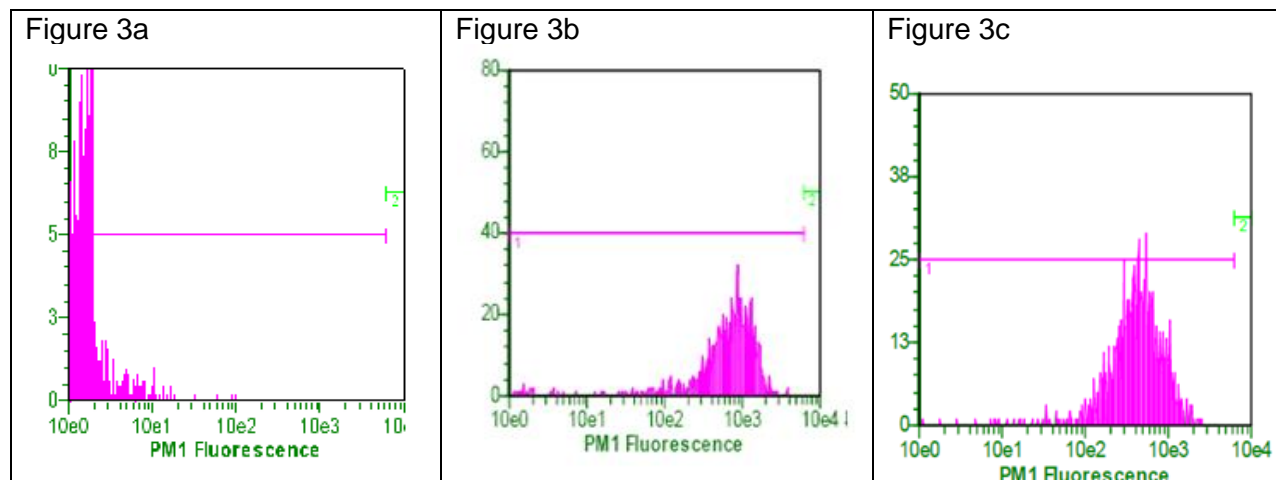
USE.

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 3. Cells staining: CFTR expression profile



CFTR expression on cell surface measured by flow cytometry (FACS) using intracellular cells staining by Streptavidin-PE.

Figure 3a: Host cells, Figure 3b: CFTR expressing cells, Figure 3c: CFTR-GFP expressing cells

References:

1. Gadsby, D.C., *et al. Nature* **440** (7083): 477–483 (2006).
2. Hillier, L.W., *et al. Nature* **424**:157-164 (2003).
3. McCann, C. M., *et al. BioTechniques* **38** (6): 945–952 (2005).

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