



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### Nav1.7 – HEK 293 Cell line

Catalog #: 60507

#### Description

Stable recombinant HEK293 cell line expressing human Nav1.7 (Genbank # Q15858) fused at the C-terminus to Green Fluorescent Protein [Ex. ~395 nm, 475 nm; em ~510 nm] and with C-terminal Streptavidin-Binding Peptide (SBP) fusion. Nav1.7 is a tetrodotoxin-sensitive voltage-gated sodium channel type IX subunit alpha (SCN9A).

#### Background

Nav1.7 is a voltage-gated sodium ion channel that in humans is encoded by the SCN9A gene. It is usually expressed at high levels in two types of neurons, the nociceptive neurons at dorsal root ganglion and trigeminal ganglion, and sympathetic ganglion neurons, which are part of the autonomic (involuntary) nervous system.

Nav1.7 is present at the endings of pain-sensing nerves, the nociceptors, close to the region where the impulse is initiated. The Nav1.7 channel produces a rapidly activating and inactivating current which is sensitive to the level of tetrodotoxin. Knockout mice that lack Nav1.7 in nociceptors showed reduced response to inflammatory pain [1].

#### Sequence

A synthetic codon-optimized DNA sequence encoding human Nav1.7 protein [2] with C-terminal Green Fluorescent Protein (GFP) and C-terminal Streptavidin-Binding Peptide (SBP) [3] tag is stably integrated in tetracycline-inducible HEK293 cells.

#### Sequence of the Nav1.7 Tag:

```
EGGGGSMVSKGEELFTGVVPIILVELDGDVNGHKFSVSGEGEGDATYGKLTLLKFICTTGKLPVPWPTLVTT  
LTYGVQCFSRYPDHMKQHDFFKSAMPEGYVQERTIFFKDDGNYKTRAEVKFEGDTLVNRIELKGIIDFKED  
GNILGHKLEYNYNVSHNVYIMADKQKNGIKVNFKIRHNIEDGSVQLADHYQQNTPIGDGPVLLPDPNHYLST  
QSALSKDPNEKRDHMLLEFVTAAGITLGMDELYKVEGLVPRGSGSLVPRGSSAKETAATAAKFERQHMDSG  
ATETSQVAPAGAAAMDEKTTGWRGGHVVEGLAGELEQLRARLEHHPQGQREP
```

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Applications

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

## Format

Each vial contains  $2 \times 10^6$  cells in 1 ml of 10% DMSO

## Host cell

HEK293 cells, tetracycline-inducible

## Recommended Storage

Immediately upon receipt, store in liquid nitrogen.

## MW

The calculated molecular weight is 242 kDa.

## Stability

The cell line has been demonstrated to be stable for at least seven continuous passages. For optimal results, it is recommended to use the cells prior to the 7<sup>th</sup> passage. Upon receipt, amplify the cells in culture and make several frozen aliquots for future use.

## Propagation Medium and Culture Conditions

**Thaw Medium 4 (BPS Cat. #60181):** DMEM/F12 (1:1) (Hyclone # SH30271.01) supplemented with 10% FBS (Life Technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone # SV30010)

**Growth Medium 4A (BPS Cat. #79535):** Thaw Medium 4 (BPS Cat. #60181) plus 10 µg/ml Blastidicin (Life Technologies # R210-01), and 200 µg/ml Zeocin (invivogen # ant-zn-1p).

Cells should be grown at 37° with 5% CO<sub>2</sub> using Growth Medium 4A (Thaw Medium 4, Blastidicin, and Zeocin).

**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 4 (**no Blastidicin and Zeocin**), spin down the cells, and resuspend cells in pre-warmed Thaw Medium 4 (**no Blastidicin and Zeocin**). Transfer resuspended cells to a T25 flask and culture in 37°C CO<sub>2</sub>

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)

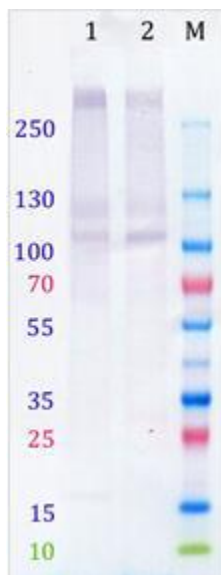
incubator. At first passage switch to Growth Medium 4A (**contains Blasticidin and Zeocin**). Cells should be split before they reach complete confluence.

**To passage the cells**, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add Growth Medium 4A and transfer to a centrifuge tube. Spin down cells, resuspend cells in Growth Medium 4A and seed appropriate aliquots of cell suspension into new culture vessels.

### Induction of the target protein expression

To express Nav1.7, cells are induced with DMEM/F12 50/50 supplemented with 10% FBS, 1% Penicillin Streptomycin, 1  $\mu$ g/ml Doxycycline (Biochemika #44577) and 3 mM Na butyrate (Acros Organics #263190250) for 24 hours prior to cell harvesting or assay.

**Figure 1. Western Blot of the Nav1.7 expressing cells.** Western Blot of HEK293 Nav1.7 cells (lanes 1, 2) stained with anti-Sodium channel Nav1.7, clone N68/6 (Millipore, Cat.No. MABN41) with follow staining with Alkaline Phosphatase conjugated Anti-mouse IgG (Rockland Immunochemicals, Cat.No. 610-1502). M: molecular weight marker.



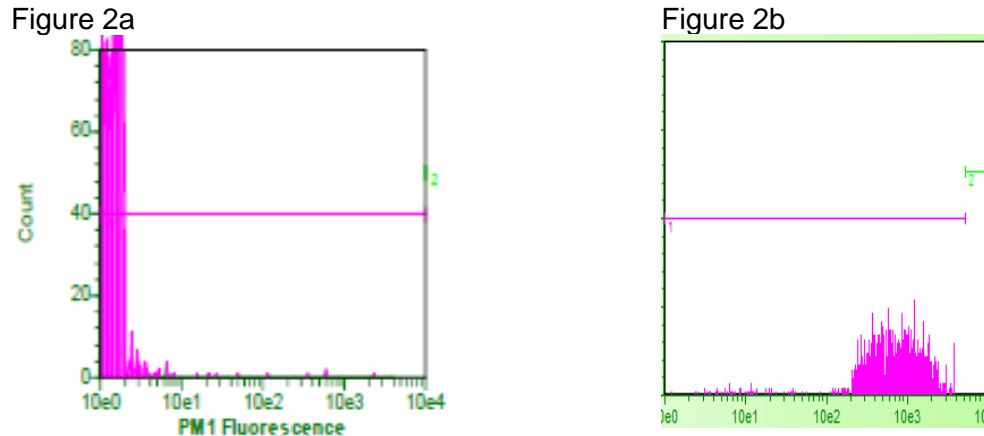
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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 2. Cell Expression Profile



NaV1.7 expression measured by flow cytometry (FACS) using intracellular staining by saponin-permeabilized cells and PE-labeled streptavidin.

Figure 2a: HEK293 Host cells, Figure 2b: NaV1.7 expressing HEK293 cells (#60507) stained with anti-NaV 1.7 clone 68/6 monoclonal antibody and PE-labeled anti-mouse.

## References

1. Catterall, W.A. *Cell and Developmental Biology* **16**: 521–555 (2000).
2. Choi, J.S., *et al. Neurology* **67**:1563-1567 (2006).
3. Li, Y., *et al. Protein Science* **20** (1): 140–149 (2011).

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

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