



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**

### **Nav1.8/ $\beta$ 2 – HEK293 Cell Line** **Catalog #: 60521** **Lot #: 160923**

#### **Description**

Stable recombinant HEK293 cell line expressing tetracycline-inducible human Nav1.8 (Genbank #NP\_006505.2) fused to Green Fluorescent Protein (GFP) [Ex. ~395 nm, 475 nm; em ~510 nm] and the human sodium channel beta 2 subunit (Genbank #NM\_004588; SCN2B). Nav1.8 is also known as tetrodotoxin-resistant voltage-gated sodium channel type X (SCN10A).

#### **Background**

Nav1.8 is a voltage gated type X, alpha subunit sodium channel which in humans is encoded by the SCN10A gene. It is expressed in nociceptors and has been proposed as a target for the development of new analgesics. Mice deficient in Nav1.8 have deficits in sensing inflammatory pain (initiated by tissue damage/inflammation) and visceral pain (initiated by damage or injury to internal organs) but not neuropathic pain. Native sodium channels are complexes composed of the pore-forming  $\alpha$  subunit and an auxiliary  $\beta$  subunit. The sodium channel beta 2 subunit, encoded by the SCN2B gene, increases current amplitude and surface expression of sodium ion channels.

#### **Sequence**

A synthetic codon-optimized DNA sequence encoding human Nav1.8 (SCN10A gene) with C-terminal GFP-tag and C-terminal Streptavidin-Binding Peptide (SBP)-tag, and the SCN2B gene encoding human  $\beta$ 2 subunit with C-terminal FLAG-tag are stably integrated in tetracycline-inducible HEK293 cells.

#### **Applications**

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

#### **Format**

Each vial contains  $\sim 1 \times 10^6$  cells in 10% DMSO solution

#### **Host cell**

HEK293 cells, tetracycline-inducible

#### **Mycoplasma testing**

The cell line is confirmed for absence of *Mycoplasma* species using the PCR-based VenorGeM<sup>®</sup> Mycoplasma Detection kit (Sigma Aldrich).

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)

### **Recommended Storage**

Immediately upon receipt, store in liquid nitrogen.

### **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 4B (BPS Cat. #79554) :** Thaw Medium 4 (BPS Cat. #60181) plus 10 µg/ml Blastidicin (Life Technologies # R210-01), 200 µg /ml Zeocin (Invivogen # ant-zn-1p), and 100 µg/ml Hygromycin B (Hyclone #SV30070.01)

Cells should be grown at 37°C with 5% CO<sub>2</sub> using complete growth medium (Thaw Medium 4, Blastidicin, Zeocin, and Hygromycin B)

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, Zeocin, and Hygromycin**), spin down the cells, and resuspend cells in pre-warmed Thaw Medium 4 (**no Blastidicin, Zeocin, and Hygromycin**). Transfer resuspended cells to a T25 flask and culture in 37°C CO<sub>2</sub> incubator. At first passage, switch to Growth Medium 4B (**contains Blastidicin, Zeocin and Hygromycin B**). 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 complete growth medium and transfer to a centrifuge tube. Spin down cells, resuspend cells in complete growth medium and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ratio: 1:20, twice per week.

### **Induction of the target protein expression**

Induce cells in DMEM/F12 (1:1), 10% FBS, 1% Penicillin Streptomycin, 1 µg/ml Doxycycline (MP Biomedicals #0219504401) and 3 mM Na butyrate (Acros Organics #263190250) for 24 hours prior to cell harvesting or assay. Cells must be induced with doxycycline and Na-butyrate for ~24 hr to express the Nav1.8.

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

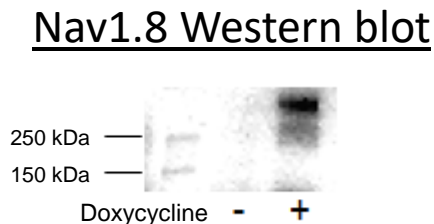
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

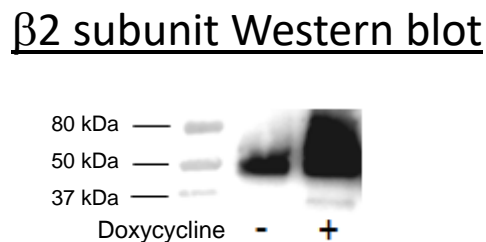
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Figure 1. Western Blot of the Nav1.8 and  $\beta$ 2 expressing the Nav1.8/ $\beta$ 2-HEK293 stable cell line.**

A) Western blot of tetracycline-induced Nav1.8-GFP, probed with anti-GFP antibody.



B) Western for  $\beta$ 2 subunit-FLAG (constitutively expressed), probed with anti-FLAG antibody.



**References**

1. Catterall, W.A. *et al. Pharmacol. Rev.* **57** (4): 397–409 (2005).
2. Muzny, D.M., *et al. Nature* **440**:1194-1198 (2006).
3. Wilson, D.S., *et al. Protein Expression and Purification* **23** (3): 440–446 (2001).

**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

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)

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)