

## PrecisION<sup>®</sup> hNav1.8/β1 Recombinant Stable Cell Line

**Catalog Number** CYL3025

**Lot Number**

See Vial

**Contents** 2 Vials, 2 x 10<sup>6</sup> to 4 x 10<sup>6</sup> in 1 mL

### Background Information

hNav1.8 is a voltage-gated sodium channel alpha subunit. It is expressed in mammalian adult dorsal root ganglion (DRG) neurons, predominantly those with small diameter cell bodies that give rise to the unmyelinated (C-type) fibres. These fibres form synapses in the dorsal horn of the spinal cord and are mainly involved in pain signaling. Additional information can be found on page 2.

### Product Information

**Description** Recombinant HEK 293 cell line expressing the human Nav1.8 (tetrodotoxin-resistant voltage-gated sodium channel type X, SCN10A) and the human sodium channel beta 1 subunit (SCN1B)

**Family** Sodium, Voltage-Gated

**Target** Nav1.8/β1

	Target Protein	Accession Number
1	Nav1.8	AF117907
2	β1	NM_001037
3	N/A	N/A
4	N/A	N/A

**Species** Human

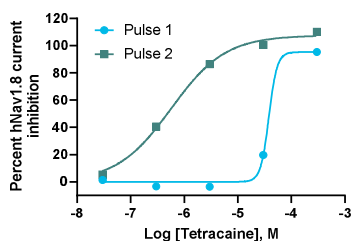
**Host Cell Type** HEK 293

**Application** Electrophysiology assay (conventional and automated patch clamp platforms)

**Storage** Vials are to be stored in vapor phase of liquid nitrogen

### Functional Performance

HEK293 cells expressing hNav1.8/β1 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



**Electrophysiology Method** QPatch

**Reference Agonist**

**Reference Antagonist** Tetracaine

**Antagonist IC<sub>50</sub> (μM)** 0.57

### Passage Stability

This cell line has been confirmed to be stable through at least 12 passages with no significant drop in assay window or change in pharmacology.

### Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

### Notes

Additional functional (pharmacological and electrophysiological) validation on multiple platforms is available upon request.

### Additional Ligand Information

**Control Compound** Tetracaine

**Vendor Name :** Sigma-Aldrich

**Vendor Catalog No.** T7383

### Additional Background Information

This channel conducts the characteristic slow, tetrodotoxin-resistant (TTX-r) currents involved in action potential initiation and transmission in these neurons and is therefore a potential target for analgesic agents. Altered expression of Nav1.8 is observed in pain models and human pain states - channel protein can be seen to re-localise from cell bodies to the site of insult or injury. There is also good evidence from antisense oligonucleotide studies for a role in development of hyperexcitable states in pain models. However KO mice show surprisingly limited pain phenotypes, possibly due to compensatory up-regulation of other Nav subtypes. Native sodium channels are multi-subunit complexes, composed of not only the pore forming  $\alpha$  subunit, but also auxiliary  $\beta$  subunits. The  $\beta$ 1 subunit is known to increase peak current amplitude as well as increasing sodium ion channel expression at the cell surface (Vijayaragavan et al., 2004).

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)

**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

**General product information:** [www.discoverx.com](http://www.discoverx.com)

### Limited Use License Agreement

**These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.**

**Please visit [discoverx.com/license](http://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.**