

PrecisION[®] hNav1.6 Recombinant Stable Cell Line

Catalog Number CYL3010

Lot Number

See Vial

Contents 2 Vials, 2 x 10⁶ to 4 x 10⁶ in 1 mL

Background Information

hNav1.6 is primarily expressed in the output neurones of the cerebellum, cortex and hippocampus of the CNS and in the Nodes of Ranvier of sensory and motor axons of the PNS. hNav1.6 is involved with the initiation and transmission of action potentials in central neurons and their myelinated axons. Mutation or disruption of the Nav1.6 gene causes cerebellar ataxia and motor end plate disease in mice. This channel is a potential target for analgesic and antiepileptic drugs.

Product Information

Description Recombinant HEK 293 cell line expressing the human Nav1.6 (type VIII voltage-gated sodium channel alpha subunit)

Family Sodium, Voltage-Gated

Target Nav1.6

	Target Protein	Accession Number
1	Nav1.6	AF050736
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

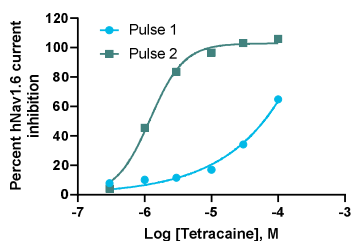
Host Cell Type HEK 293

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 hNav1.6 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



Electrophysiology Method QPatch

Reference Agonist ATX-II

Reference Antagonist Tetracaine

Antagonist IC₅₀ (μM) 1.22

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

N/A

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com

Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com

General product information: 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 for a list of products that are governed by limited use label license terms and relevant patent and trademark information.