

PrecisION[®] hHCN2 Recombinant Stable Cell Line

Catalog Number CYL3041

Lot Number

See Vial

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

Background Information

Hyperpolarization-activated, cyclic nucleotide-modulated (HCN) channels contribute to rhythmic spontaneous electrical activity in the heart and CNS. Interference with HCN2 activity can lead to various forms of cardiac arrhythmias, due to dysfunction of the sinus node. The predominant role of HCN2 in these cells may be the stabilization of a normal diastolic membrane potential.

Product Information

Description Recombinant HEK 293 cell line expressing the human hyperpolarization-activated, cyclic nucleotide-gated potassium channel 2 (HCN2)

Family HCN, Hyperpolarized

Target HCN2

	Target Protein	Accession Number
1	HCN2	NM_001194
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

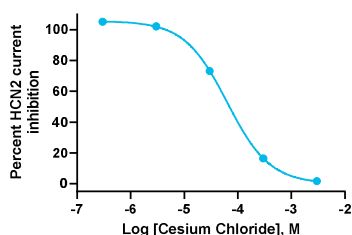
Host Cell Type HEK 293

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 hHCN2 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



Electrophysiology Method QPatch

Reference Agonist

Reference Antagonist Cesium Chloride

Antagonist IC₅₀ (μM) 63.50

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 Cesium Chloride

Vendor Name : Tocris

Vendor Catalog No. 4739

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.