

PrecisION[®] hKv3.3 Recombinant Stable Cell Line

Catalog Number CYL3045

Lot Number

See Vial

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

Background Information

Kv3.3 are voltage-gated, fast inactivating Shaw channels which play an important role in regulating neuronal output by modulating the delay and frequency of high frequency neuronal firing. They are highly expressed in areas of the brain involved with motor control such as the cerebellum. They have also been shown to be expressed at the mammalian neuromuscular junction where they are thought to regulate transmitter release.

Product Information

Description Recombinant CHO-K1 cell line expressing the human voltage-gated potassium channel Kv3.3

Family Potassium, Voltage-Gated

Target Kv3.3

	Target Protein	Accession Number
1	Kv3.3	NM_004977
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

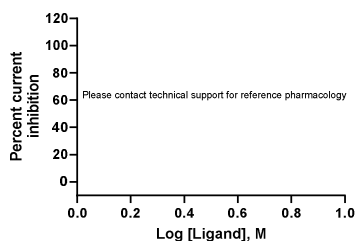
Host Cell Type CHO-K1

Application Electrophysiology assay (conventional and automated patch clamp platforms)

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

Functional Performance

CHO-K1 cells expressing hKv3.3 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



Electrophysiology Method MPC

Reference Agonist

Reference Antagonist 4-AP

Antagonist IC₅₀ (μM)

Passage Stability

Please contact technical support.

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 4-AP

Vendor Name : Sigma-Aldrich

Vendor Catalog No. 275875

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.