

PrecisION[®] hKv1.7 Recombinant Stable Cell Line

Catalog Number CYL3020

Lot Number

See Vial

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

Background Information

Kv1.7 is a voltage-gated, delayed rectifier, potassium channel. It is expressed in the heart, skeletal muscle and pulmonary arteries. As this channel has properties similar to the ultra-rapidly activating (IKur) current of the heart, it is potentially a candidate channel for involvement in human cardiac disorders.

Product Information

Description Recombinant CHO-K1 cell line expressing the human voltage-gated potassium channel Kv1.7

Family Potassium, Voltage-Gated

Target Kv1.7

	Target Protein	Accession Number
1	Kv1.7	NM_031886
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

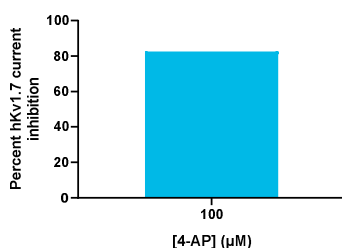
Host Cell Type CHO-K1

Application Electrophysiology assay (conventional and automated patch clamp platforms)

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

Functional Performance

CHO cells expressing hKv1.7 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

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 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.