

## PrecisION<sup>®</sup> hKv1.3 Recombinant Stable Cell Line

**Catalog Number** CYL3016

**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

Kv1.3 is widely expressed throughout the CNS and in organs/cells involved with immune system such as spleen, tonsils, T-lymphocytes, B-lymphocytes and macrophages. Kv1.3 is involved in the regulation of the resting membrane potential and calcium signalling in lymphocytes and oligodendrocytes. This ion channel is a target for immunosuppressants and Kv1.3 inhibitors have been proven effective treatment in slowing the progression of multiple sclerosis.

### Product Information

**Description** Recombinant CHO-K1 cell line expressing the human Kv1.3 potassium channel

**Family** Potassium, Voltage-Gated

**Target** Kv1.3

	Target Protein	Accession Number
1	Kv1.3	NM_002232
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

**Species** Human

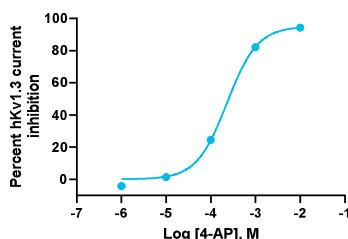
**Host Cell Type** CHO-K1

**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 hKv1.3 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



**Electrophysiology Method** QPatch

**Reference Agonist**

**Reference Antagonist** 4-AP

**Antagonist IC<sub>50</sub> (μM)** 231.50

### 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** 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](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.**