

## PrecisION<sup>®</sup> hERG Recombinant Stable Cell Line

**Catalog Number** CYL3038

**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

The FDA, ICH S7B guidelines set out a non-clinical testing strategy to assess the effects of pharmaceuticals on ventricular repolarisation and proarrhythmic risk. The most common mechanism for this is inhibition of the delayed rectified potassium current IKr. The human ether-à-go-go-related gene (hERG) encodes the potassium ion channel responsible for IKr.

### Product Information

**Description** Recombinant CHO-K1 cell line expressing the human ERG (ether-a-go-go related gene) potassium channel

**Family** Potassium, Voltage-Gated

**Target** hERG

	Target Protein	Accession Number
1	hERG	U04270
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

**Species** Human

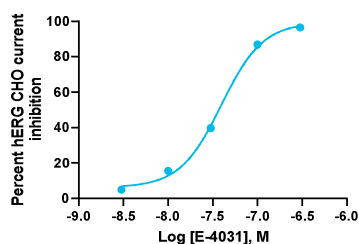
**Host Cell Type** CHO-K1

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



**Electrophysiology Method** QPatch

**Reference Agonist**

**Reference Antagonist** E-4031

**Antagonist IC<sub>50</sub> (μM)** 0.04

### 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** E-4031

**Vendor Name :** Sigma-Aldrich

**Vendor Catalog No.** M5060

### 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.**