

## PrecisION<sup>®</sup> hCav3.2 Recombinant Stable Cell Line

**Catalog Number** CYL3075

**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 T-type calcium channel (hCav3.2) is a type of voltage-dependent calcium channel, the  $\alpha 1$  subunit determines most of the channel's properties. The T-type calcium channel is one of the components that produce the pacemaker potential in the SA node of the heart. T-type calcium channel blockers are used primarily as antiepileptics. Additional information can be found on page 2.

### Product Information

**Description** Recombinant HEK 293 cell line expressing the human T-type calcium channel Cav3.2

**Family** Calcium, Voltage-Gated

**Target** Cav3.2

	Target Protein	Accession Number
1	Cav3.2	NM_021098
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

**Species** Human

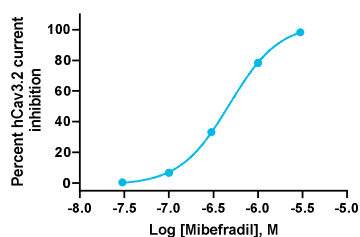
**Host Cell Type** HEK 293

**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 hCav3.2 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



**Electrophysiology Method** QPatch

**Reference Agonist**

**Reference Antagonist** Mibefradil

**Antagonist IC<sub>50</sub> (μM)** 4.80

### 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** Mibefradil

**Vendor Name :** Tocris

**Vendor Catalog No.** 2198

### Additional Background Information

Two voltage-dependent calcium channels play critical roles in the physiology of cardiac muscle: L-type calcium channel ('L' for Long-lasting) and T-type calcium channels ('T' for Transient) voltage-gated calcium channels. L-type channels are important in sustaining an action potential, while T-type channels are important in initiating them.

**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.**