

## PrecisION<sup>®</sup> hGABAA $\alpha$ 5/ $\beta$ 3/ $\gamma$ 2 Recombinant Stable Cell Line

**Catalog Number** CYL3073

**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

Gamma-aminobutyric acid (GABA)-gated ion channels are widely distributed in the mammalian brain and are major mediators of inhibitory synaptic transmission. A typical GABA<sub>A</sub> ion channel has a pentameric structure consisting of 5 protein subunits, often  $\alpha$ ,  $\beta$  and  $\gamma$ , combining to form a central ion conducting pore across the cell membrane. Additional information can be found on page 2.

### Product Information

**Description** Recombinant HEK 293 cell line expressing the human GABAA  $\alpha$ 5,  $\beta$ 3 and  $\gamma$ 2 ion channel subunits

**Family** Chloride, Ligand-Gated

**Target** GABAA  $\alpha$ 5/ $\beta$ 3/ $\gamma$ 2

	Target Protein	Accession Number
1	GABAA $\alpha$ 5	NM_000810
2	GABAA $\beta$ 3	NM_000814
3	GABAA $\gamma$ 2	NM_000816
4	N/A	N/A

**Species** Human

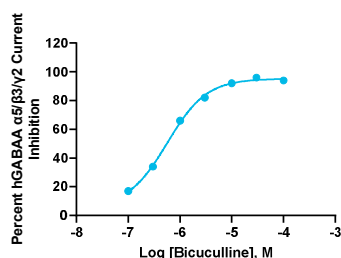
**Host Cell Type** HEK 293

**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 hGABAA  $\alpha$ 5/ $\beta$ 3/ $\gamma$ 2 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



**Electrophysiology Method** IonFlux

**Reference Agonist** GABAA

**Reference Antagonist** Bicuculline

**Antagonist IC<sub>50</sub> ( $\mu$ M)** 0.57

### 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** Bicuculline

**Vendor Name :** Tocris

**Vendor Catalog No.** 0130

### Additional Background Information

In humans there are six genes that encode  $\alpha$  subunits, three that encode  $\beta$ , three that encode  $\gamma$ , and an additional seven genes that encode other subunits whose function is less-well understood than the  $\alpha$ ,  $\beta$  and  $\gamma$  subunits. GABA<sub>A</sub> ion channels open and close in response to secretion of GABA from presynaptic terminals. GABA<sub>A</sub>  $\alpha 5$  channels are typically comprised of  $\alpha 5$ ,  $\beta 3$ , and  $\gamma 2$  subunits and this combination is preferentially expressed in the mammalian hippocampus (Wisden et al., 1992). Much pharmaceutical research interest has focused on  $\alpha 5$ -selective "inverse agonists" at the benzodiazepine (BZ) site, which decrease the current elicited by a concentration of GABA that produces a 20% response, i.e. an EC<sub>20</sub> (Atack et al. 2006, Dawson et al. 2006). These compounds show promise as dementia treatments because in animal models they enhance cognition with minimal proconvulsant or anxiogenic side-effects.

**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.