

## PrecisION<sup>®</sup> hnAChR $\alpha$ 3/ $\beta$ 4 Recombinant Stable Cell Line

**Catalog Number** CYL3057

**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

hnAChR $\alpha$ 3 $\beta$ 4 receptor is a nicotinic acetylcholine receptor located in the autonomic ganglia and adrenal medulla parts of the brain. Its activation causes excitation in the pre-and post-synapse.

### Product Information

**Description** Recombinant HEK 293 cell line expressing the  $\alpha$ 3 and  $\beta$ 4 subunits of the human nicotinic acetylcholine receptor

**Family** Ligand-Gated

**Target** nAChR  $\alpha$ 3/ $\beta$ 4

	Target Protein	Accession Number
1	nAChR $\alpha$ 3	M37981
2	$\beta$ 4	NM_000750
3	N/A	N/A
4	N/A	N/A

**Species** Human

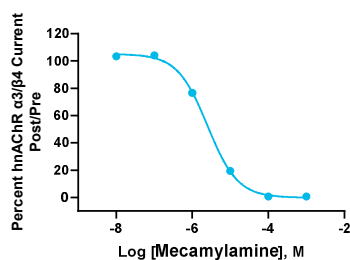
**Host Cell Type** HEK 293

**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 hnAChR  $\alpha$ 3/ $\beta$ 4 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



**Electrophysiology Method** IonFlux

**Reference Agonist** Acetylcholine

**Reference Antagonist** Mecamylamine

**Antagonist IC<sub>50</sub> ( $\mu$ M)** 2.52

### 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** Mecamylamine

**Vendor Name :** Tocris

**Vendor Catalog No.** 2843

### 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.