

PrecisION[®] hnAChR α 1/ β 1/ δ / ϵ Recombinant Stable Cell Line

Catalog Number CYL3052

Lot Number

See Vial

Contents 2 Vials, 2 x 10⁶ to 4 x 10⁶ in 1 mL

Background Information

The α 1/ β 1/ δ / ϵ receptor is the major nicotinic receptor found in adult mammalian muscle. These muscle nACh receptors are involved in the fast synaptic transmission at the skeletal-muscular junction and their activation causes muscular contraction. This cell line would be useful for safety counter screening compounds for any undesirable action on this nicotinic receptor subtype.

Product Information

Description Recombinant HEK 293 cell line expressing the four subunits α 1, β 1, δ and ϵ of the human nicotinic acetylcholine receptor type I

Family Ligand-Gated

Target nAChR α 1/ β 1/ δ / ϵ

	Target Protein	Accession Number
1	nAChR α 1	NM_000079
2	nAChR β 1	NM_000747
3	nAChR δ	NM_000751
4	nAChR ϵ	NM_000080

Species Human

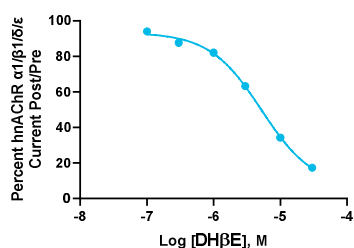
Host Cell Type HEK 293

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 α 1/ β 1/ δ / ϵ were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



Electrophysiology Method IonFlux

Reference Agonist Acetylcholine

Reference Antagonist DHβE

Antagonist IC₅₀ (μM) 5.21

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 DH β E

Vendor Name : Tocris

Vendor Catalog No. 2349

Additional Background Information

N/A

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com

Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com

General product information: 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 for a list of products that are governed by limited use label license terms and relevant patent and trademark information.