

PrecisION[®] hGluR6 Recombinant Stable Cell Line

Catalog Number CYL3049

Lot Number

See Vial

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

Background Information

hGluR6 is a member of the ligand-gated, kainate family of glutamate receptors also known as GRIK2. It can form both hetero and homo-tetrameric, cation permeable, channels and is expressed at high levels in the hippocampus and striatum of the CNS. Pre- and postsynaptic kainate receptors are able to regulate both neurotransmitter release and excitability in a synapse-specific manner. Additional information can be found on page 2.

Product Information

Description Recombinant HEK 293 cell line expressing the human GluR6 ion channel

Family Ligand-Gated

Target GluR6

	Target Protein	Accession Number
1	GluR6	NM_021956
2	N/A	N/A
3	N/A	N/A
4	N/A	N/A

Species Human

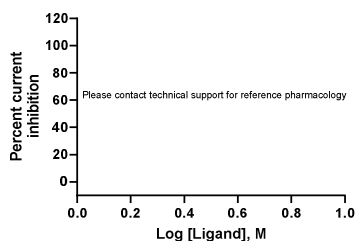
Host Cell Type HEK 293

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 hGluR6 were characterized in terms of their pharmacological and biophysical properties using whole-cell patch clamp techniques.



Electrophysiology Method

Reference Agonist

Reference Antagonist

Antagonist IC₅₀ (μM)

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

Vendor Name :

Vendor Catalog No.

Additional Background Information

GRIK2 has been implicated in a range of neurophysiological conditions such as autism, obsessive compulsive disorder, as well as disease states e.g. Schizophrenia, epilepsy and Huntingdon's disease.

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.