

cAMP Hunter™ eXpress GRM6 CHO-K1 GPCR Assay

Catalog Number: 95-0105E2

Lot Number: See Vial

Contents: 3.75 x 10⁶ cells per vial in 0.2 mL

Background

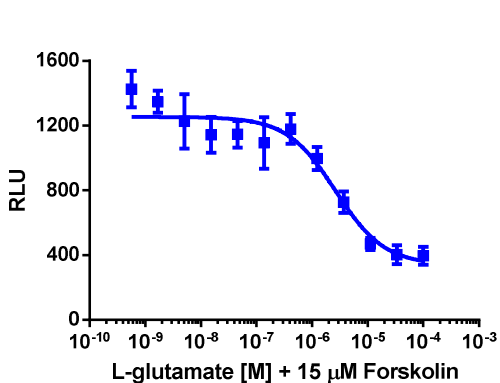
cAMP Hunter eXpress cells are engineered to overexpress naturally Gi or Gs coupled, wildtype GPCRs and are designed to detect changes in intracellular cAMP levels in response to agonist stimulation of the receptor. These cells have been modified to prevent long term propagation and expansion using a proprietary compound that has no apparent effect on assay performance.

Product Information

Target GPCR:	GRM6
Description:	Metabotropic glutamate receptor 6
Receptor Family:	Metabotropic glutamate
Coupling:	Gi/Go
Accession Number:	NM_000843
GPCR Species:	Human
Cell Type:	CHO-K1
Storage:	Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

Functional Performance

Cells were plated in a 96-well plate and stimulated with a control agonist (+ Forskolin for Gi targets), using the assay conditions described below. Following stimulation, signal was detected according to the recommended protocol. Please refer below for information on control compounds.



Cell Number/Well:	30000
Control Agonist:	L-glutamate
Cell Plating Reagent:	AssayComplete™ Cell Plating 24 Reagent
Cell Incubation Time (Hours):	24
Agonist Incubation Time (minutes):	30
Agonist Incubation Temperature (°C):	37
EC₅₀ for Agonist Stimulation (nM):	2644
Signal:Background at Agonist E_{max}:	3.5

Important! This assay requires an additional step: After cell plating reagent is replaced with buffer (3 reagent addition protocol) or cAMP antibody working solution (2 reagent addition protocol), pre-incubate cells for 1 hour at 37° C, prior to addition of ligand. Then continue according to recommended protocol.

Additional Ligand Information

Control Agonist: L-glutamate

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1197)

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.

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