

cAMP Hunter™ eXpress PROKR1 CHO-K1 GPCR Assay

Catalog Number: 95-0090E2

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: PROKR1

Description: Prokineticin receptor 1

Receptor Family: Prokineticin

Coupling: Gs & Gq

Accession Number: NM_138964

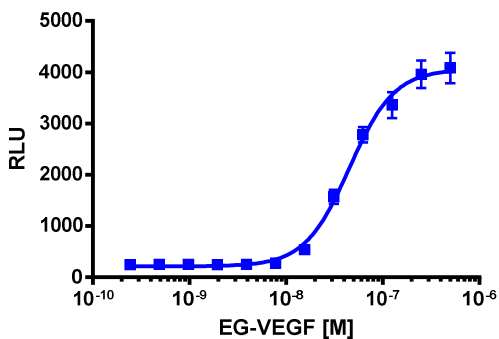
GPCR Species: Human

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: Recombinant Human EG-VEGF

Cell Plating Reagent: AssayComplete™ Cell Plating 2 Reagent

Cell Incubation Time (Hours): 24

Agonist Incubation Time (minutes): 30

Agonist Incubation Temperature (°C): 37

EC₅₀ for Agonist Stimulation (nM): 45

Signal:Background at Agonist E_{max}: 18.9

Additional Ligand Information

Control Agonist: Recombinant Human EG-VEGF

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1022)

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