

PathHunter® eXpress CHRM4 CHO-K1 β -Arrestin GPCR Assay

Catalog Number: 93-0349E2

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

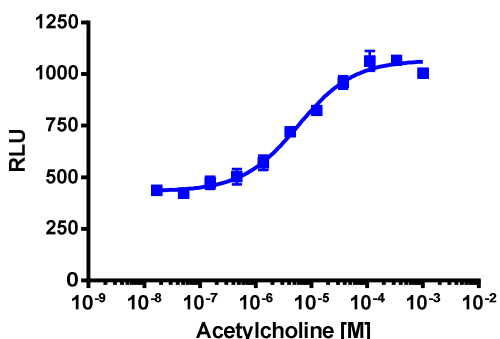
PathHunter eXpress β -Arrestin GPCR cells are engineered to co-express the ProLink™ (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged β -Arrestin. Activation of the GPCR-PK induces β -Arrestin-EA recruitment, forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal. 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: | CHRM4 |
| Description: | Cholinergic receptor, muscarinic 4 |
| Receptor Family: | Acetylcholine |
| Coupling: | Gs & Gi/Go |
| Accession Number: | NM_000741.2 |
| GPCR Species: | Human |
| β-Arrestin Isoform: | β -Arrestin-2 |
| ProLink™ Tag: | PK1 |
| 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, 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: | 10000 |
| Control Agonist: | Acetylcholine |
| Cell Plating Reagent: | AssayComplete™ Cell Plating 0 Reagent |
| Cell Incubation Time (Hours): | 24 |
| Agonist Incubation Time (Minutes): | 90 |
| Agonist Incubation Temperature (°C): | 37 |
| EC₅₀ for Agonist Stimulation (nM): | 5743 |
| Signal:Background at Agonist E_{max}: | 2.5 |

Additional Ligand Information

Control Agonist: Acetylcholine

Vendor: Eurofins DiscoverX® (Catalog No. 92-1125)

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