

PathHunter® eXpress GCGR Total GPCR Internalization Assay

Catalog Number: 93-0820E1 **Lot Number:** See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

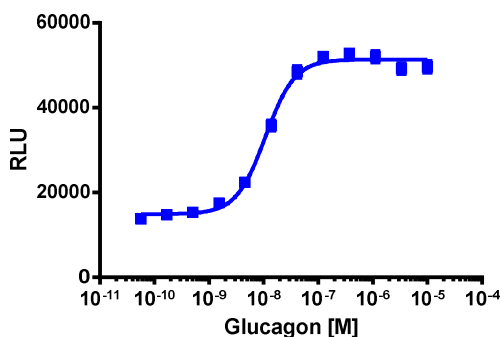
PathHunter eXpress Total GPCR Internalization cells are engineered to co-express the Enzyme Acceptor (EA) tagged GPCR, and a ProLink™ (PK) tag localized to the endosomes. Activation of the GPCR-EA induces internalization of the receptor in PK-tagged endosomes, forcing complementation of the two β-galactosidase 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: GCGR
Description: Glucagon receptor
Receptor Family: Glucagon
Coupling: Gs
Accession Number: NM_000160
GPCR Species: Human
Cell Type: HEK 293
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: Glucagon
Cell Plating Reagent: AssayComplete™ Cell Plating 0 Reagent
Cell Incubation Time (Hours): 24
Agonist Incubation Time (minutes): 180
Agonist Incubation Temperature (°C): 37
EC₅₀ for Agonist Stimulation (nM): 10.8
Signal:Background at Agonist E_{max}: 3.6

Additional Ligand Information

Control Agonist: Glucagon

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1087)

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