

PathHunter[®] eXpress GLP2R CHO-K1 β -Arrestin GPCR Assay

Catalog Number: 93-0572E2

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

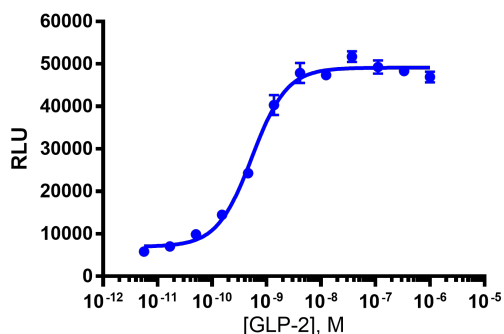
PathHunter eXpress β -Arrestin GPCR cells are engineered to co-express the ProLink™ (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged β -Arrestin. Activation of the GPCR-PK induces β -Arrestin-EA recruitment, forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The resulting 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:	GLP2R
Description:	Glucagon-like peptide receptor 2
Receptor Family:	Glucagon
Coupling:	Gs
Accession Number:	NM_004246.1
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:	Glucagon-like Peptide 2 (GLP-2)
Cell Plating Reagent:	AssayComplete™ Cell Plating 2 Reagent
Cell Incubation Time (Hours):	24
Agonist Incubation Time (Minutes):	90
Agonist Incubation Temperature (°C):	37
EC₅₀ for Agonist Stimulation (nM):	0.54
Signal:Background at Agonist E_{max}:	8.1

Additional Ligand Information

Control Agonist: Glucagon-like Peptide 2 (GLP-2)

Vendor: Eurofins DiscoverX® (Catalog No. 92-1079)

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