

## PathHunter® eXpress GLP1R Activated GPCR Internalization Assay

**Catalog Number:** 93-0724E3

**Lot Number:**

See Vial

**Contents:** 1 x 10<sup>6</sup> cells per vial in 0.1 mL

### Background

PathHunter eXpress Activated GPCR Internalization cells are engineered to co-express an untagged GPCR, Enzyme Acceptor (EA) tagged  $\beta$ -Arrestin, and a ProLink™ (PK) tag localized to the endosomes. Activation of the GPCR induces  $\beta$ -Arrestin recruitment, followed by internalization of the Receptor/Arrestin-EA complex in PK-tagged endosomes. This forces complementation of the two  $\beta$ -galactosidase enzyme fragments (EA and PK), forming functional enzyme that 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:** GLP1R

**Description:** Glucagon-like peptide receptor 1

**Receptor Family:** Glucagon

**Coupling:** Gs

**Accession Number:** NM\_002062.3

**GPCR Species:** Human

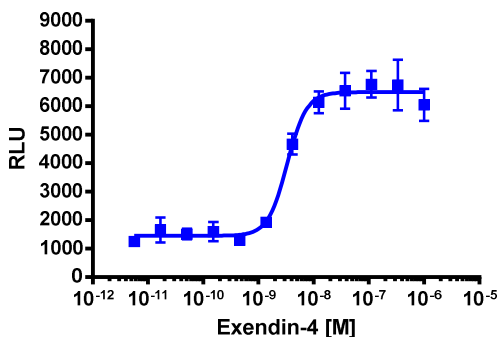
**$\beta$ -Arrestin Isoform:**  $\beta$ -Arrestin-2

**Cell Type:** U2OS

**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:**

Exendin-4

**Cell Plating Reagent:**

AssayComplete™ Cell Plating 0 Reagent

**Cell Incubation Time (Hours):**

**Agonist Incubation Time (minutes):**

180

**Agonist Incubation Temperature (°C):**

37

**EC<sub>50</sub> for Agonist Stimulation (nM):**

3.3

**Signal:Background at Agonist E<sub>max</sub>:**

4.8

**Important!** This assay requires an additional step: Please refer to Additional Protocol Information section.

### Additional Ligand Information

**Control Agonist:** Exendin-4

**Vendor:** Eurofins DiscoverX<sup>®</sup> (Catalog No. 92-1115)

### Additional Protocol Information

It is recommended that the control compound be serially diluted in a base media (or buffer) + 1% FBS to ensure optimal assay performance.

### 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](https://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](mailto:CustomerServiceDRX@eurofins.com)

**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

**General product information:** [www.discoverx.com](https://www.discoverx.com)