

## 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

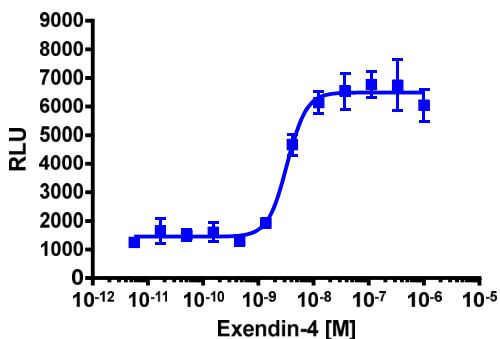
PathHunter eXpress Activated GPCR Internalization cells are engineered to co-express an untagged GPCR, Enzyme Acceptor (EA) tagged β-Arrestin, and a ProLink™ (PK) tag localized to the endosomes. Activation of the GPCR induces β-Arrestin recruitment, followed by internalization of the Receptor/Arrestin-EA complex in PK-tagged endosomes. This forces complementation of the two β-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  
**β-Arrestin Isoform:** β-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):** 48  
**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® (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 <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.