

## PathHunter® eXpress GLP1R Total GPCR Internalization Assay

**Catalog Number:** 93-1075E3

**Lot Number:**

See Vial

**Contents:** 1.2 x 10<sup>6</sup> cells per vial in 0.1 mL

### Background

PathHunter eXpress Total GPCR Internalization cells are engineered to co-express the ProLink™ (PK) tagged GPCR, and an Enzyme Acceptor (EA) tag localized to the endosomes. Activation of the GPCR-PK induces internalization of the receptor in EA-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:** GLP1R

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

**Receptor Family:** Glucagon

**Coupling:** Gs

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

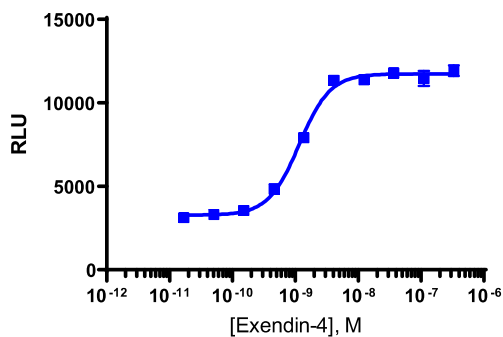
**GPCR Species:** Human

**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 5 Reagent

**Cell Incubation Time (Hours):** 48

**Agonist Incubation Time (minutes):** 180

**Agonist Incubation Temperature (°C):** 37

**EC<sub>50</sub> for Agonist Stimulation (nM):** 1.1

**Signal:Background at Agonist E<sub>max</sub>:** 3.6

### Additional Ligand Information

**Control Agonist:** Exendin-4

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1115)

### 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.