

## PathHunter® eXpress GHSR U2OS $\beta$ -Arrestin-1 GPCR Assay

**Catalog Number:** 93-0242E3

**Lot Number:** See Vial

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

### Background

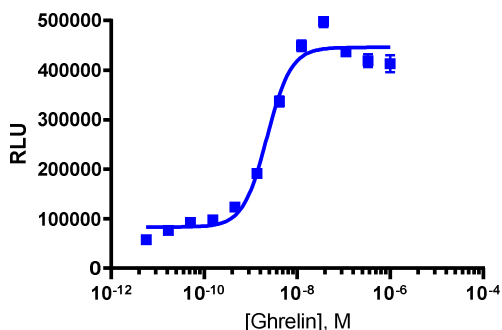
PathHunter eXpress  $\beta$ -Arrestin GPCR cells are engineered to co-express the ProLink™ (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged  $\beta$ -Arrestin. Activation of the GPCR-PK induces  $\beta$ -Arrestin-EA recruitment, forcing complementation of the two  $\beta$ -galactosidase enzyme fragments (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

<b>Target GPCR:</b>	GHSR
<b>Description:</b>	Ghrelin receptor
<b>Receptor Family:</b>	Ghrelin
<b>Coupling:</b>	Gq
<b>Accession Number:</b>	NM_198407
<b>GPCR Species:</b>	Human
<b><math>\beta</math>-Arrestin Isoform:</b>	$\beta$ -Arrestin-1
<b>ProLink™ Tag:</b>	PK1
<b>Cell Type:</b>	U2OS
<b>Storage:</b>	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.



<b>Cell Number/Well:</b>	10000
<b>Control Agonist:</b>	Ghrelin
<b>Cell Plating Reagent:</b>	AssayComplete™ Cell Plating 5 Reagent
<b>Cell Incubation Time (Hours):</b>	48
<b>Agonist Incubation Time (Minutes):</b>	90
<b>Agonist Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for Agonist Stimulation (nM):</b>	2.3
<b>Signal:Background at Agonist E<sub>max</sub>:</b>	7.6

### Additional Ligand Information

**Control Agonist:** Ghrelin

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1345)

### 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)