

## PathHunter® CHO-K1 BDKRB2 $\beta$ -Arrestin Cell Line

**Catalog Number:** 93-0189C2 **Lot Number:** See Vial  
**Contents:** 2 vials, 1 x 10<sup>6</sup> cells per vial in 1 mL

### Background

PathHunter  $\beta$ -Arrestin GPCR cell lines 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 enzyme fragments. The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal. PathHunter cell lines expressing Gq-coupled receptors can also be used to detect calcium mobilization.

### Product Information

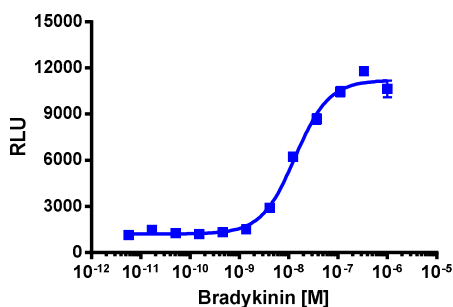
**Target GPCR:** BDKRB2  
**Description:** Bradykinin receptor b2  
**Receptor Family:** Bradykinin  **$\beta$ -Arrestin Isoform:**  $\beta$ -Arrestin-2  
**Accession Number:** NM\_000623.3 **ProLink™ Tag:** PK1  
**Coupling:** Gq **Cell Type:** CHO-K1  
**GPCR Species:** Human  
**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

### Functional Performance

Cells were plated in a 384-well plate and incubated overnight at 37°C and 5% CO<sub>2</sub> to allow the cells to attach and grow. Cells were then stimulated with a control agonist using the assay conditions described below. Following stimulation, signal from Arrestin recruitment was detected using the PathHunter Detection Kit according to the recommended protocol. Calcium mobilization was detected using the Calcium No Wash<sup>PLUS</sup> detection kit. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.

#### PathHunter Arrestin Assay

**Cell Number/Well:** 5000  
**Agonist Incubation Time (minutes):** 90  
**Agonist Incubation Temp. (°C):** 37



**EC<sub>50</sub> for Arrestin Recruitment (nM):** 13.5

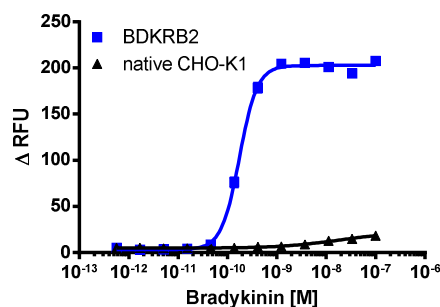
**Signal:Background at agonist E<sub>max</sub>:** 10.3

**Control Antagonist:** HOE 140 **Arrestin Assay IC<sub>50</sub> (nM):** 61.9

**Note:** This cell line was developed and quality control tested via the PathHunter Arrestin Assay only. Calcium mobilization assay was run independently from the PathHunter Arrestin Assay. Calcium data is background subtracted and represented as  $\Delta$ RFU (Relative Fluorescence Units).

#### Ca<sup>2+</sup> Mobilization Assay

**Cell Number/Well:** 10000  
**Signal Read Time (@ 2 sec intervals):** 2 minutes  
**Signal Read Temperature (°C):** RT



**EC<sub>50</sub> for Calcium Mobilization (nM):** 0.17

### Passage Stability

This cell line has been confirmed to be stable through 10 passages with no significant drop in assay window or change in EC<sub>50</sub>. Passage stability testing was conducted using the PathHunter Arrestin Assay only.

### Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

### Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	PathHunter® Detection Kit	93-0001
Ca <sup>2+</sup> Detection	Calcium No Wash <sup>PLUS</sup>	90-0091
Cell Culture	AssayComplete™ Cell Culture Kit-107	92-3107G
Cell Plating	AssayComplete™ Cell Plating 2 Reagent	93-0563R2A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T2	92-4102TR
Cell Freezing	AssayComplete™ Freezing Reagent F2	92-5102FR

\*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

### Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	300	92-0029
AssayComplete™ G418	800	92-0030

### Additional Ligand Information

**Control Compound:** Bradykinin

**Vendor:** DiscoverX® (Catalog No. 92-1053)

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail [info@discoverx.com](mailto:info@discoverx.com). For additional information, please visit [discoverx.com](http://discoverx.com).

## 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.