

# PathHunter<sup>®</sup> CHO-K1 MRGPRE $\beta$ -Arrestin Orphan GPCR Cell Line

**Catalog Number:** 93-0438C2A **Lot Number:** See Vial  
**Contents:** 2 vials, 1 x 10<sup>6</sup> cells per vial in 1 mL

## Background

PathHunter  $\beta$ -Arrestin Orphan 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  $\beta$ -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

## Product Information

**Target GPCR:** MRGPRE  
**Description:** MAS-related GPR, member E  
**Receptor Family:** Class A Orphan  
**Accession Number:** NM\_001039165.2  
**GPCR Species:** Human  
 **$\beta$ -Arrestin Isoform:**  $\beta$ -Arrestin-2  
**ProLink™ Tag:** PK1  
**Cell Type:** CHO-K1  
**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

## Functional Performance

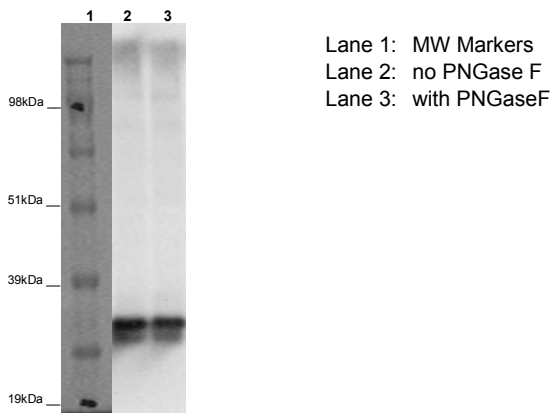


Figure 1. Cell lysates prepared from PathHunter  $\beta$ -Arrestin Orphan GPCR cell lines were treated with PNGase F (Glyko; Cat. #GKE-5003), run on a SDS-PAGE gel and analyzed. Untreated lane resolves a band of appropriate size corresponding to GPCR-PK fusion protein and the PNGase F treated lane resolves a deglycosylated band indicative of proper expression and folding of GPCR protein.

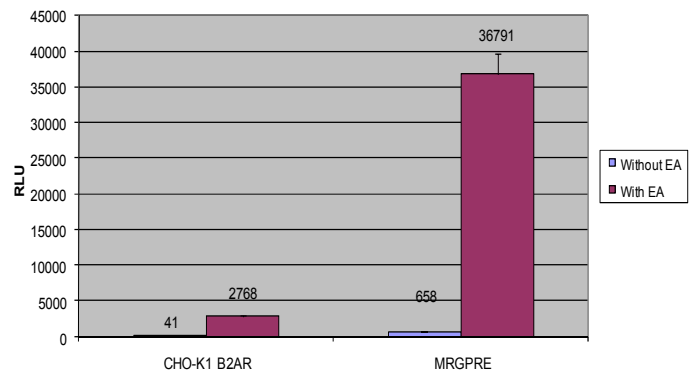


Figure 2. PathHunter  $\beta$ -Arrestin Orphan GPCR cells were analyzed for basal activity as well as GPCR-ProLink™ expression by comparing the ratio of signal between untreated cells and cells treated with saturating amounts of exogenous EA, using ProLink™ Detection Kit (DrX: 92-0006). Signal from complementation of ProLink™ and EA fragments correlates to the amount of GPCR-PK expression in the cell line.

### Passage Stability

This cell line has been confirmed to be stable through 10 passages with no significant change in GPCR-PK expression level.

### 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 <sup>®</sup> Detection Kit  | 93-0001        |
| Cell Culture    | AssayComplete™ Cell Culture Kit-107    | 92-3107G       |
| Cell Plating    | AssayComplete™ Cell Plating 1 Reagent  | 93-0563R1A     |
| 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        |

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.