

# PathHunter® eXpress TRHR CHO-K1 $\beta$ -Arrestin GPCR Assay

**Catalog Number:** 93-0239E2

**Lot Number:** See Vial

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

## Background

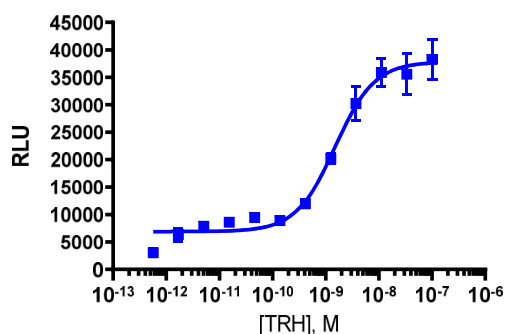
PathHunter eXpress  $\beta$ -Arrestin GPCR cells are engineered to co-express the ProLink™ (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged  $\beta$ -Arrestin. Activation of the GPCR-PK induces  $\beta$ -Arrestin-EA recruitment, forcing complementation of the two  $\beta$ -galactosidase enzyme fragments (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>                         | TRHR  |
| <b>Description:</b>                         | Thyrotropin-releasing hormone receptor  |
| <b>Receptor Family:</b>                     | Thyrotropin releasing hormone   |
| <b>Coupling:</b>                            | Gq  |
| <b>Accession Number:</b>                    | AY493373  |
| <b>GPCR Species:</b>                        | Human   |
| <b><math>\beta</math>-Arrestin Isoform:</b> | $\beta$ -Arrestin-2   |
| <b>ProLink™ Tag:</b>                        | PK1   |
| <b>Cell Type:</b>                           | CHO-K1  |
| <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>                              | TRH                                   |
| <b>Cell Plating Reagent:</b>                         | AssayComplete™ Cell Plating 0 Reagent |
| <b>Cell Incubation Time (Hours):</b>                 | 24                                    |
| <b>Agonist Incubation Time (Minutes):</b>            | 90                                    |
| <b>Agonist Incubation Temperature (°C):</b>          | 37                                    |
| <b>EC<sub>50</sub> for Agonist Stimulation (nM):</b> | 1.5                                   |
| <b>Signal:Background at Agonist E<sub>max</sub>:</b> | 12.5                                  |

### Additional Ligand Information

**Control Agonist:** TRH

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1072)

### 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)