

## PathHunter<sup>®</sup> eXpress cINSRb Functional Assay

**Catalog Number:** 93-1018E3

**Lot Number:**

See Vial

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

### Background

PathHunter RTK Functional cell lines are engineered to co-express a ProLink™ (PK) tagged RTK, an Enzyme Acceptor (EA) tagged SH2 domain, and, in some cases, an untagged co-receptor. Activation of the RTK-PK induces receptor dimerization, leading to SH2-EA recruitment, and 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 RTK:** cINSRb

**RTK Accession #:** XM\_005587740.1

**Description:** Insulin receptor, isoform B

**Kinase Species:** Macaca fascicularis

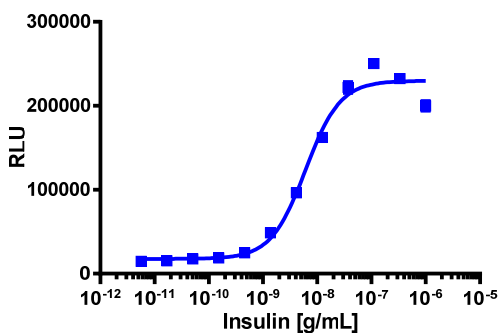
**SH2 Domain:** PLCG1

**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:**

Insulin, Human

**Cell Plating Reagent:**

AssayComplete™ Cell Plating 16 Reagent

**Cell Incubation Time (Hours):**

48

**Agonist Incubation Time (minutes):**

180

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

RT

**EC<sub>50</sub> for Agonist Stimulation (ng/mL):**

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

### Additional Information

**Control Agonist:** Insulin, Human

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1088)

### 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](http://www.discoverx.com)