

# PathHunter® eXpress rhTrkA-rhP75 Functional Assay

**Catalog Number:** 93-1009E3

**Lot Number:**

See Vial

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

## Background

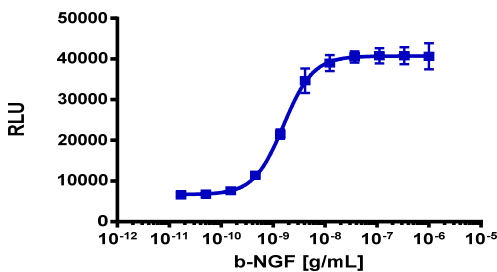
PathHunter RTK Functional cell lines are engineered to co-express a ProLink™ (PK) tagged RTK, an Enzyme Acceptor (EA) tagged SH2 domain, and, in some cases, an untagged co-receptor. Activation of the RTK-PK induces receptor dimerization, leading to SH2-EA recruitment, and forcing complementation of the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

<b>Target RTK:</b>	rhTrkA	<b>Co-receptor:</b>	rhP75
<b>RTK Accession #:</b>	XM_001114243.2	<b>Accession #:</b>	Customer sequence
<b>Description:</b>	Neurotrophic tyrosine kinase receptor type 1	<b>Description:</b>	nerve growth factor receptor
<b>Kinase Species:</b>	Rhesus		
<b>SH2 Domain:</b>	SHC1		
<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.



**Cell Number/Well:**

10000

**Control Agonist:**

Recombinant Human b-NGF

**Cell Plating Reagent:**

AssayComplete™ Cell Plating 16 Reagent

**Cell Incubation Time (Hours):**

24

**Agonist Incubation Time (minutes):**

180

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

RT

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

1.6

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

6.2

### Additional Information

**Control Agonist:** Recombinant Human b-NGF

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1023)

### 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)