

PathHunter[®] eXpress IGF1R Functional Assay

Catalog Number: 93-0505E1 **Lot Number:** See Vial
Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

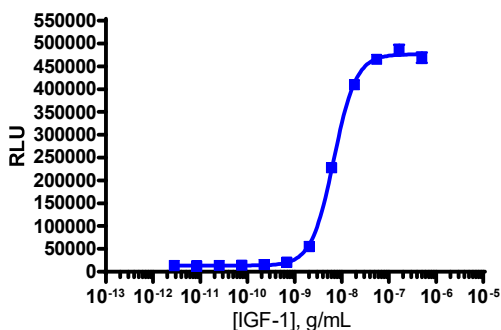
PathHunter eXpress Total RTK Internalization cells are engineered to co-express the ProLink™ (PK) tagged RTK, and an Enzyme Acceptor (EA) tag localized to the endosomes. Activation of the RTK-PK induces internalization of the receptor in EA-tagged endosomes, forcing complementation of the two β-galactosidase 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: IGF1R
Description: Insulin-Like Growth Factor 1 Receptor
Accession Number: NM_000875
RTK Species: Human
Cell Type: HEK 293
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: Recombinant Human IGF-1
Cell Plating Reagent: AssayComplete™ Cell Plating 17 Reagent
Cell Incubation Time (Hours): 0
Agonist Incubation Time (minutes): 180
Agonist Incubation Temperature (°C): RT
EC₅₀ for Agonist Stimulation ng/mL 6.8
Signal:Background at Agonist E_{max}: 35.1

Additional Ligand Information

Control Agonist: Recombinant Human IGF-1
Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1030)

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 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
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com