

PathHunter® eXpress c-Ret-GFRα1 Functional Assay

Catalog Number: 93-1054E3

Lot Number:

See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

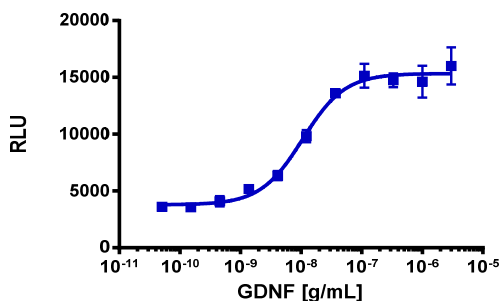
PathHunter RTK Functional cell lines are engineered to co-express a ProLink™ (PK) tagged RTK, an Enzyme Acceptor (EA) tagged SH2 domain, and, in some cases, an untagged co-receptor. Activation of the RTK-PK induces receptor dimerization, leading to SH2-EA recruitment, and forcing complementation of the 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:	c-Ret	Co-receptor:	GFRα1
RTK Accession #:	NM_020975.4	Accession #:	NM_001145453.1
Description:	Proto-oncogene tyrosine-protein kinase receptor Ret	Description:	GDNF family receptor alpha-1
Kinase Species:	Human		
SH2 Domain:	SHC1		
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:	GDNF
Cell Plating Reagent:	AssayComplete™ Cell Plating 16 Reagent
Cell Incubation Time (Hours):	4
Agonist Incubation Time (minutes):	960
Agonist Incubation Temperature (°C):	37
EC₅₀ for Agonist Stimulation (ng/mL):	10.6
Signal:Background at Agonist E_{max}:	4.4

Additional Information

Control Agonist: GDNF

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1299)

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