

Certificate Of Analysis

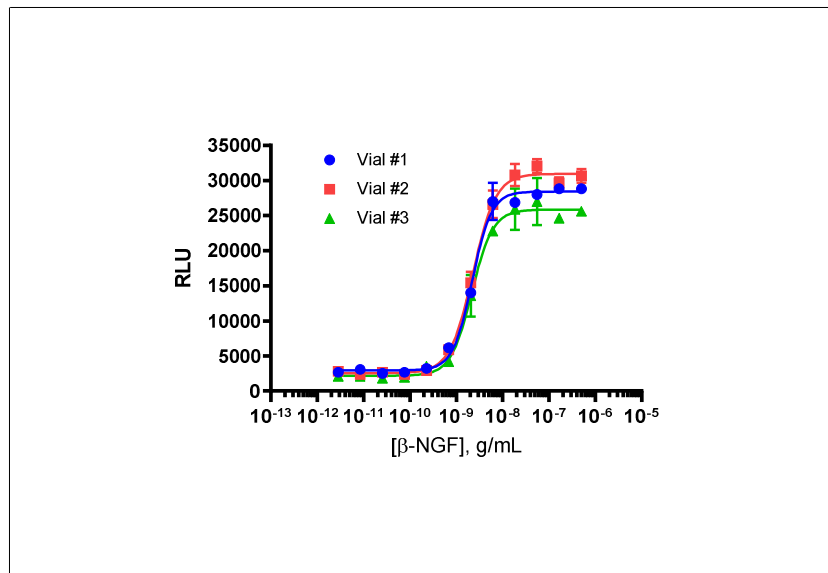
Background

PathHunter[®] RTK Bioassay Cells 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 receptor phosphorylation, and 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.

Product Name	PathHunter[®] U2OS TrkA-P75 Bioassay Cells
Cryovial Label	U2OS TrkA-P75 Bioassay Cells
Bioassay Catalog #	93-0529Y3
Bioassay Manufactured Lot #	20C2404
Passage # @ Freezing	4

Assay Information	
Target	TrkA
Target Accession Number	NM_002529
Target Description	Neurotrophic tyrosine kinase receptor type 1
Target 2	P75
Target 2 Accession Number	NM_002507
Target 2 Description	nerve growth factor receptor
SH2 Domain	SHC1
Species	Human
Cell	U2OS
CP Reagent	AssayComplete [™] Cell Plating 16 Reagent (DiscoverX, 93-0563R16A)
Ligand	Recombinant Human b-NGF (DiscoverX, 92-1023)
Ligand Diluent	AssayComplete [™] Protein Dilution Buffer
Detection Kit	PathHunter [®] Detection Kit (DiscoverX, 93-0001M)
Cell Number/Well	5,000
Cell Seeding Time (hours)	24
Ligand Inc Time (minutes)	180
Agonist Inc Temperature (°C)	RT

Cell Density Information	
Cell Number (millions)	1.2
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	96
Recovery After 48 Hours (%)	174
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 10.8
	Vial 2 11.0
	Vial 3 12.2
EC ₅₀ (g/mL)	Vial 1 2.2 x 10 ⁻⁹
	Vial 2 2.3 x 10 ⁻⁹
	Vial 3 2.2 x 10 ⁻⁹



Shipping and Storage Information	
Shipping Conditions	Dry Ice
Storage Conditions	Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen.
Manufacturing Date	March 2020
Expiration Date	September 2023

Shelf life of over 3 years has been established for DiscoverX cell lines and Assay-Ready Cells in general, when stored in the vapor phase of liquid nitrogen.

Documented by / Date: _____

Approved by / Date: _____