

## Certificate Of Analysis

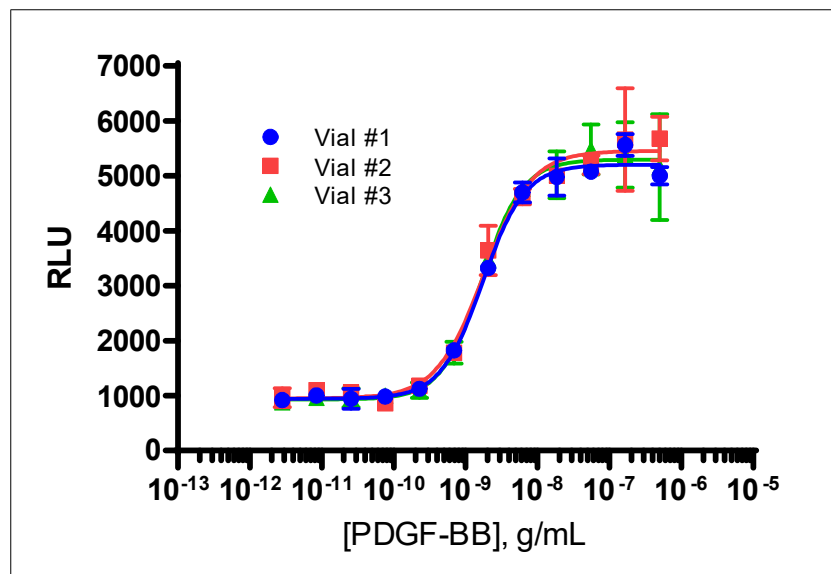
### Background

PathHunter<sup>®</sup> RTK Bioassay Cells are engineered to co-express a ProLink<sup>™</sup> (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  $\beta$ -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

<b>Product Name</b>	<b>PathHunter<sup>®</sup> U2OS PDGFRA Bioassay</b>
Cryovial Label	U2OS PDGFRA Functional Kinase Bioassay Cells
Bioassay Catalog #	93-0823Y3
Bioassay Manufactured Lot #	21K2811
Passage # @ Freezing	4

<b>Assay Information</b>	
Target	PDGFRA
Target Accession Number	NM_006206.4
Target Description	Platelet-derived growth factor receptor alpha
Target 2	N/A
Target 2 Accession Number	N/A
Target 2 Description	N/A
SH2 Domain	PLCG1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete <sup>™</sup> Cell Plating 22 Reagent (DiscoverX, 93-0563R22A)
Ligand	Recombinant Human PDGF-BB (DiscoverX, 92-1110)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter <sup>®</sup> Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	10,000
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 (%)	94
Recovery After 48 Hours (%)	211
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 5.4
	Vial 2 5.9
	Vial 3 5.9
EC <sub>50</sub> (g/mL)	Vial 1 1.7 X 10 <sup>-9</sup>
	Vial 2 1.8 X 10 <sup>-9</sup>
	Vial 3 1.7 X 10 <sup>-9</sup>



<b>Shipping and Storage Information</b>	
Shipping Conditions	Dry Ice
Storage Conditions	Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen.
Manufacturing Date	November 2021
Expiration Date	May 2025

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: \_\_\_\_\_