

## Certificate Of Analysis

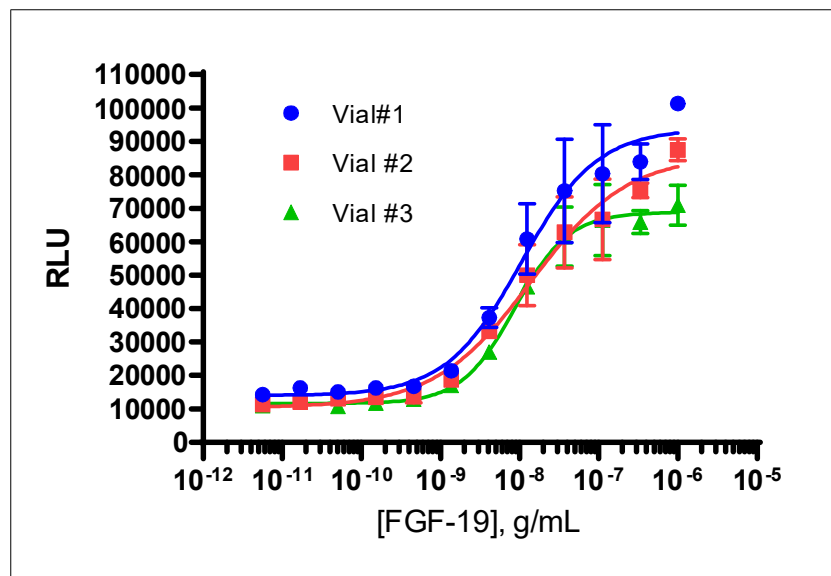
### Background

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 FGFR4-<math>\beta</math>-Klotho Bioassay Cells</b>
Cryovial Label	U2OS FGFR4-Beta-Klotho Functional Kinase Bioassay Cells
Bioassay Catalog #	93-1060Y3
Bioassay Manufactured Lot #	22B2807
Passage # @ Freezing	4

<b>Assay Information</b>	
Target	FGFR4
Target Accession Number	NM_002011
Target Description	Fibroblast growth factor receptor 4
Target 2	$\beta$ -Klotho
Target 2 Accession Number	NM_175737.3
Target 2 Description	KLB
SH2 Domain	PLCG1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete <sup>™</sup> Cell Plating 0 Reagent (DiscoverX, 93-0563R0A)
Ligand	FGF-19 (DiscoverX, 92-1301)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter <sup>®</sup> Detection Kit (DiscoverX, 93-0001)
Cell Number/Well	10,000
Cell Seeding Time (hours)	0
Ligand Inc Time (minutes)	960
Agonist Inc Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	1.2
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	97
Recovery After 48 Hours (%)	152
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 7.1
	Vial 2 7.8
	Vial 3 6.4
EC <sub>50</sub> (g/mL)	Vial 1 10.5 x 10 <sup>-9</sup>
	Vial 2 14.0 x 10 <sup>-9</sup>
	Vial 3 8.7 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	March 2022
Expiration Date	September 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: \_\_\_\_\_