

Certificate Of Analysis

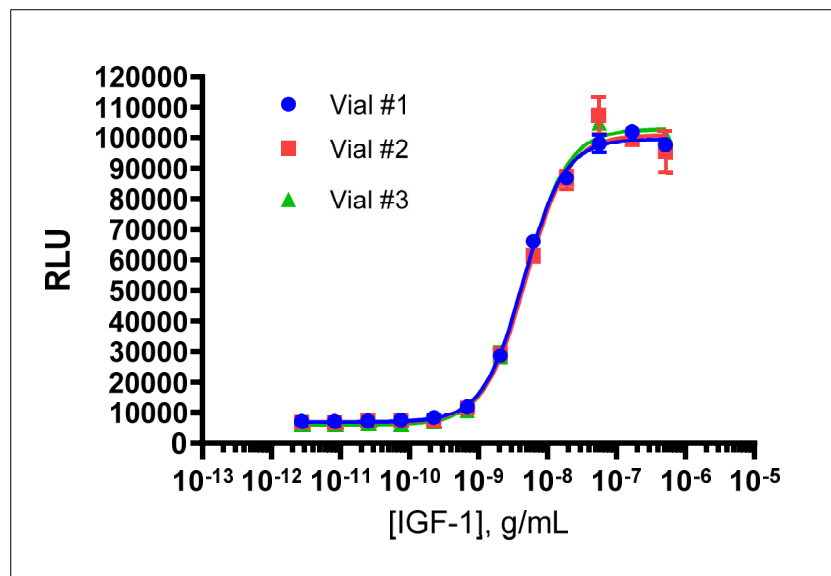
Background

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[®] HEK 293 IGF1R Bioassay
Cryovial Label	HEK 293 IGF1R Functional Kinase Bioassay Cells
Bioassay Catalog #	93-0505Y1
Bioassay Manufactured Lot #	21E2703
Passage # @ Freezing	3

Assay Information	
Target	IGF1R
Target Accession Number	NM_000875
Target Description	Insulin-like growth factor 1 receptor
Target 2	N/A
Target 2 Accession Number	N/A
Target 2 Description	N/A
SH2 Domain	SHC1
Target Species	Human
Cell	HEK 293
CP Reagent	AssayComplete [™] Cell Plating 17 Reagent (DiscoverX, 93-0563R17A)
Ligand	Recombinant Human IGF-1 (DiscoverX, 92-1030)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933)
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 (%)	94
Recovery After 48 Hours (%)	305
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 13.7
	Vial 2 13.8
	Vial 3 15.8
EC ₅₀ (g/mL)	Vial 1 4.5 x 10 ⁻⁹
	Vial 2 4.9 x 10 ⁻⁹
	Vial 3 4.8 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	June 2021
Expiration Date	December 2024

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: _____