

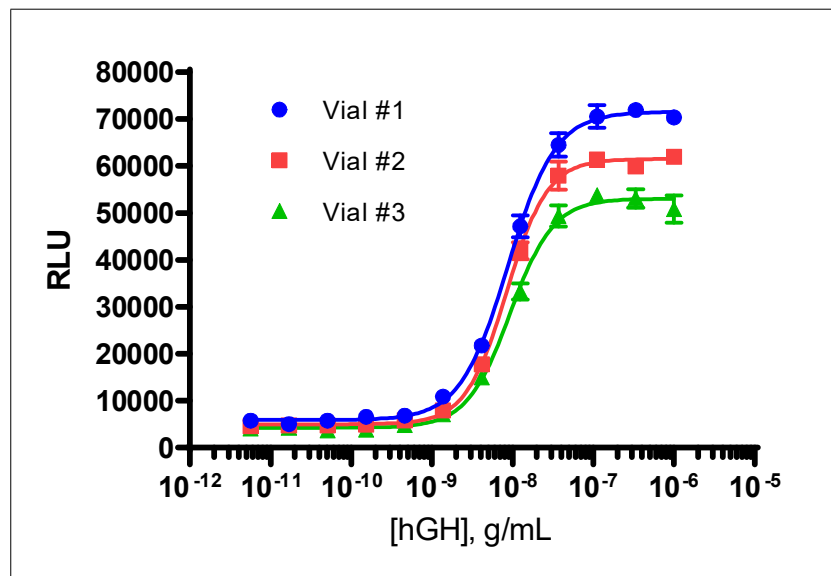
Certificate Of Analysis

Background

PathHunter[®] Cytokine Bioassay Cells are engineered to co-express a ProLink[™] (PK) tagged Cytokine receptor, an untagged CTK, and an Enzyme Acceptor (EA) tagged SH2 domain. Activation of the Receptor-PK induces phosphorylation of the receptor by the CTK. This leads to SH2-EA recruitment, 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 GHR-JAK1 Bioassay Cells
Cryovial Label	U2OS GHR-JAK1 Kinase Bioassay Cells
Bioassay Catalog #	93-0756Y3
Bioassay Manufactured Lot #	22E3109
Passage # @ Freezing	6
Assay Information	
Target 1	GHR
Target 1 Accession Number	N/A
Target 1 Description	Growth hormone receptor
Target 2	JAK1
Target 2 Accession Number	NM_002227
Target 2 Description	Janus kinase 1
SH2 Domain	PLCG1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete [™] Cell Plating 4 Reagent (DiscoverX, 93-0563R4A)
Ligand	Recombinant Human Growth Hormone (DiscoverX, 92-1338)
Ligand Diluent	DB-B5
Detection Kit	PathHunter [®] 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)	4.0
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	97
Recovery After 48 Hours (%)	173
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 12.2
	Vial 2 13.5
	Vial 3 12.5
EC ₅₀ (g/mL)	Vial 1 8.7 x 10 ⁻⁹
	Vial 2 8.4 x 10 ⁻⁹
	Vial 3 9.3 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 2022
Expiration Date	June 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: _____