

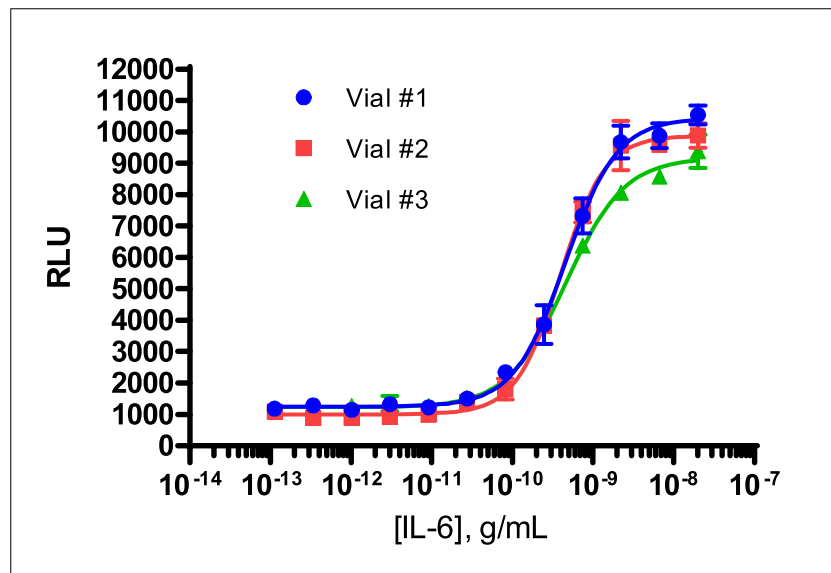
Certificate Of Analysis

Background

This PathHunter® Bioassay detects ligand induced dimerization of two subunits of a receptor-dimer pair. The cells have been engineered to co-express one receptor subunit fused to Enzyme Donor (ED), and a second dimer partner fused to Enzyme Acceptor (EA). Binding of an agonist to one receptor subunit induces it to interact with its dimer partner, forcing complementation of the two enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes a substrate to generate a chemiluminescent signal.

Product Name	PathHunter® U2OS IL6R/IL6ST Bioassay
Cryovial Label	U2OS IL6R/IL6ST Dimerization Bioassay Cells
Bioassay Catalog Number	93-1045B3
Bioassay Manufactured Lot #	22D2606
Passage # @ Freezing	4
Assay Information	
Target 1	IL6R
Target 1 Amino Acid Range	1 - 468
Target 1 Accession Number	NM_000565.3
Target 1 Description	interleukin 6 receptor
Target 1 Tag	EA
Target 2	IL6ST
Target 2 Amino Acid Range	1 - 647
Target 2 Accession Number	NM_002184.3
Target 2 Description	interleukin 6 signal transducer
Target 2 Tag	PK1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete™ Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	IL-6 (DiscoverX, 92-1285)
Ligand Diluent	AssayComplete™ Cell Plating 5 Reagent
Detection Kit	PathHunter® Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	10,000
Cell Seeding Time (hours)	24
Ligand Inc Time (minutes)	360
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 (%)	96
Recovery After 48 Hours (%)	285
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 8.9
	Vial 2 9.1
	Vial 3 7.9
EC ₅₀ (g/mL)	Vial 1 0.46 x 10 ⁻⁹
	Vial 2 0.39 x 10 ⁻⁹
	Vial 3 0.47 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	May 2022
Expiration Date	May 2026

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: _____