

Certificate Of Analysis

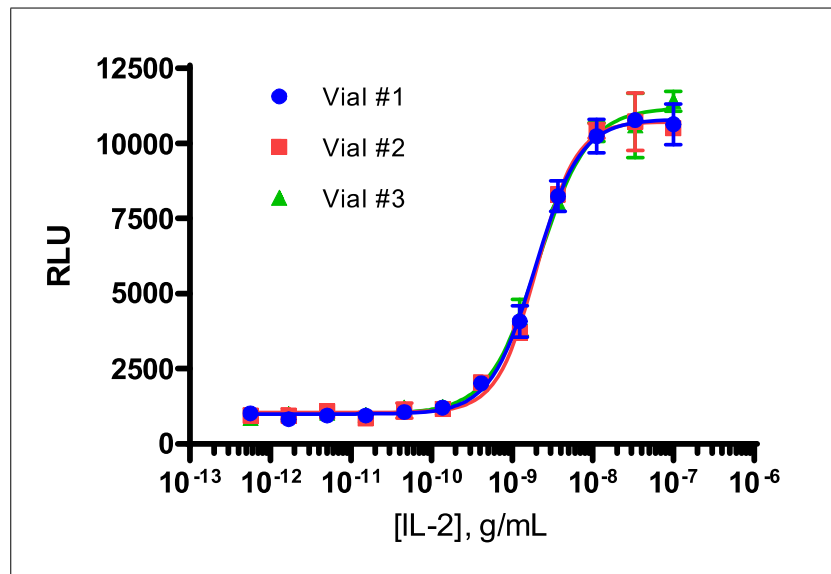
Background

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). Intracellular catalytic domains may have been deleted from the receptor subunits. 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 IL2RB/IL2RG/IL2RA Dimerization Bioassay
Cryovial Label	U2OS IL2RB/IL2RG/IL2RA Dimerization Bioassay Cells
Bioassay Catalog #	93-1003Y3
Bioassay Manufactured Lot #	22C3104
Passage # @ Freezing	4
Assay Information	
Target 1	IL2RB
Target 1 Amino Acid Range	1 - 270
Target 1 Accession Number	NM_000878.3
Target 1 Description	Interleukin 2 receptor, beta
Target 1 Tag	PK1
Target 2	IL2RG
Target 2 Amino Acid Range	1 - 290
Target 2 Accession Number	NM_000206.2
Target 2 Description	Interleukin 2 receptor, gamma
Target 2 Tag	EA
Target 3 (untagged)	IL2RA
Target 3 Amino Acid Range	1 - 272
Target 3 Accession Number	NM_000417.2
Target 3 Description	Interleukin 2 receptor, alpha
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete™ Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	IL-2 (DiscoverX, 92-1253)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	5,000
Cell Seeding Time (hours)	24
Ligand Inc Time (minutes)	360
Agonist Inc Temperature (°C)	37

Certificate Of Analysis

Cell Density Information	
Cell Number (millions)	0.6
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	96
Recovery After 48 Hours (%)	173
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 10.4
	Vial 2 11.2
	Vial 3 13.3
EC ₅₀ (g/mL)	Vial 1 1.94 x 10 ⁻⁹
	Vial 2 2.03 x 10 ⁻⁹
	Vial 3 2.11 x 10 ⁻⁹



Certificate Of Analysis

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	April 2022
Expiration Date	April 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: _____