

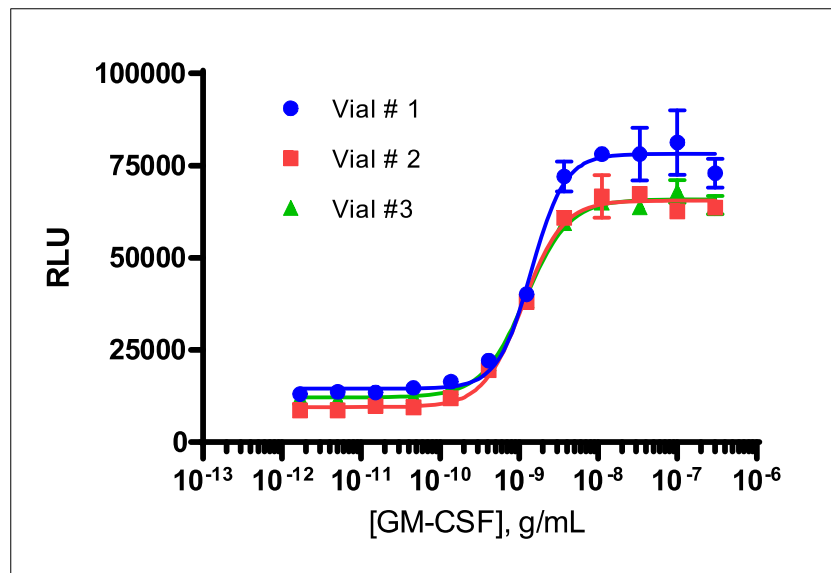
Certificate Of Analysis

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 CSF2RB/CSF2RA Bioassay Cells
Cryovial Label	U2OS CSF2RB/CSF2RA Dimerization Bioassay Cells
Bioassay Catalog Number	93-1078Y3
Bioassay Manufactured Lot #	22A0703
Passage # @ Freezing	5
Assay Information	
Target 1	CSF2RB
Target 1 Amino Acid Range	1 - 470
Target 1 Accession Number	NM_000395.2
Target 1 Description	CSF2RB protein
Target 1 Tag	PK1
Target 2	CSF2RA
Target 2 Amino Acid Range	1 - 355
Target 2 Accession Number	NM_006140.4
Target 2 Description	Granulocyte-macrophage colony-stimulating factor receptor subunit alpha
Target 2 Tag	EA
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete™ Cell Plating 22 Reagent (DiscoverX, 93-0563R22A)
Ligand	GM-CSF (DiscoverX, 92-1317)
Ligand Diluent	AssayComplete™ Dilution Buffer B5
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	5,000
Cell Seeding Time (hours)	4
Ligand Inc Time (minutes)	960
Agonist Inc Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	0.6
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	92
Recovery After 48 Hours (%)	221
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 5.6
	Vial 2 7.4
	Vial 3 5.4
EC ₅₀ (g/mL)	Vial 1 1.4 x10 ⁻⁹
	Vial 2 1.1 x10 ⁻⁹
	Vial 3 1.2 x10 ⁻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	January 2022
Expiration Date	January 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: _____