

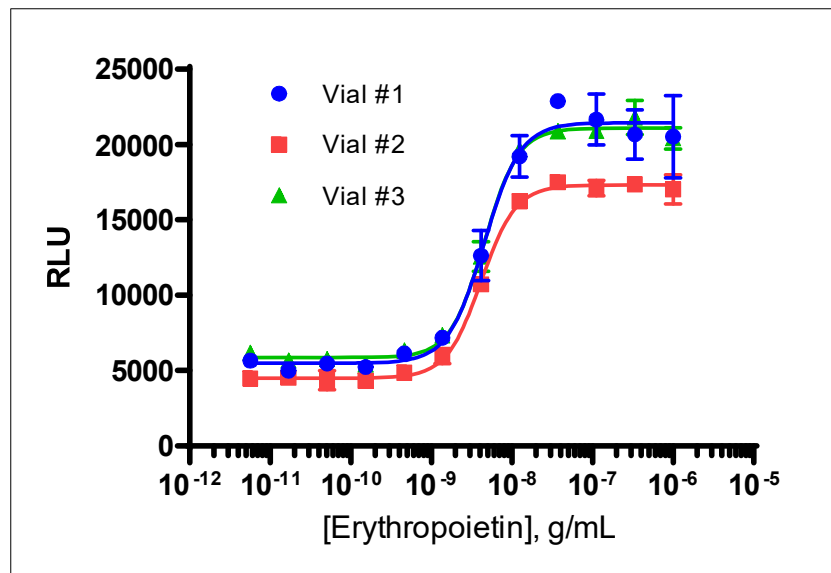
Certificate Of Analysis

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 EpoR-EpoR Bioassay
Cryovial Label	U2OS EpoR-EpoR Dimerization Bioassay Cells
Bioassay Catalog Number	93-0965Y3
Bioassay Manufactured Lot #	22J0607
Passage # @ Freezing	5
Assay Information	
Target 1	EpoR
Target 1 Amino Acid Range	1 - 508
Target 1 Accession Number	NM_000121.3
Target 1 Description	Erythropoietin receptor
Target 1 Tag	PK1
Target 2	EpoR
Target 2 Amino Acid Range	1 - 297
Target 2 Accession Number	NM_000121.3
Target 2 Description	Erythropoietin receptor
Target 2 Tag	EA
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete™ Cell Plating 0 Reagent (DiscoverX, 93-0563R0A)
Ligand	Erythropoietin (DiscoverX, 92-1224)
Ligand Diluent	Protein Dilution Buffer
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 (%)	93
Recovery After 48 Hours (%)	109
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 3.6
	Vial 2 3.8
	Vial 3 3.3
EC ₅₀ (g/mL)	Vial 1 4.5 x 10 ⁻⁹
	Vial 2 4.1 x 10 ⁻⁹
	Vial 3 4.5 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	September 2022
Expiration Date	September 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: _____