

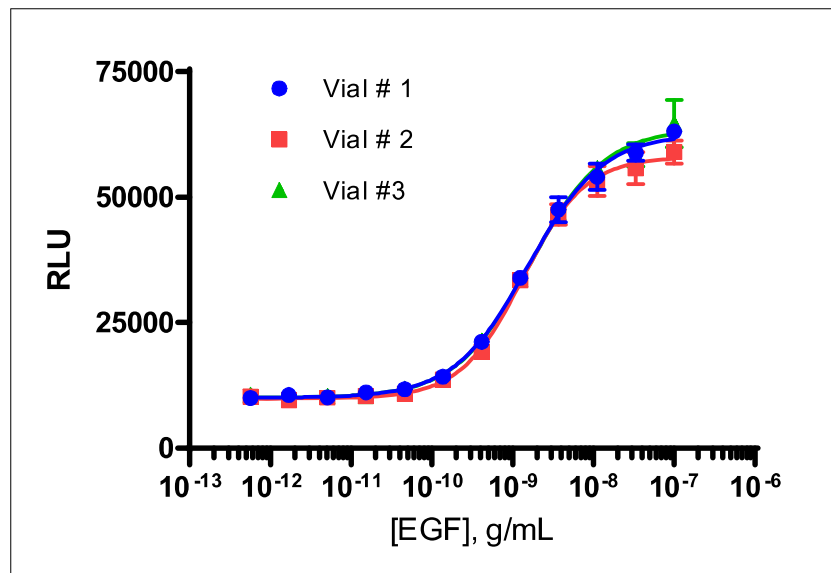
Certificate Of Analysis

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 EGFR/ErbB2 Bioassay
Cryovial Label	U2OS EGFR/ErbB2 Dimerization Bioassay Cells
Bioassay Catalog Number	93-1051Y3
Bioassay Manufactured Lot #	22C2207
Passage # @ Freezing	4
Assay Information	
Target 1	EGFR
Target 1 Amino Acid Range	1 - 679
Target 1 Accession Number	NM_005228.3
Target 1 Description	Epidermal growth factor receptor
Target 1 Tag	EA
Target 2	ErbB2
Target 2 Amino Acid Range	1 - 686
Target 2 Accession Number	NM_004448.3
Target 2 Description	erb-b2 receptor tyrosine kinase 2
Target 2 Tag	PK1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete [™] Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	Recombinant Human EGF (DiscoverX, 92-1113)
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)	1.2
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	96
Recovery After 48 Hours (%)	152
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 6.3
	Vial 2 5.8
	Vial 3 6.1
EC ₅₀ (g/mL)	Vial 1 1.6 x 10 ⁻⁹
	Vial 2 1.3 x 10 ⁻⁹
	Vial 3 1.7 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	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: _____