

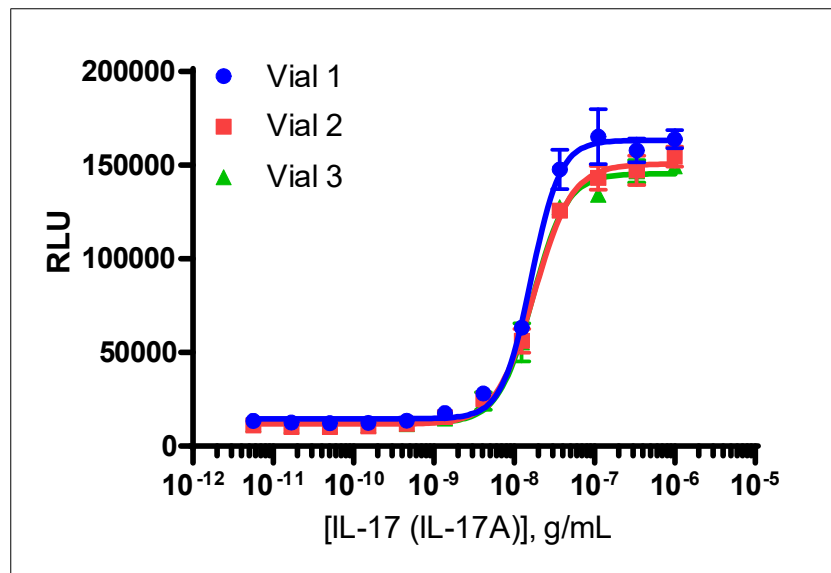
Certificate Of Analysis

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 IL17RA/IL17RC Bioassay Cells
Cryovial Label	U2OS IL17RA/IL17RC Dimerization Bioassay Cells
Bioassay Catalog Number	93-0999Y3
Bioassay Manufactured Lot #	23C0806
Passage # @ Freezing	4
Assay Information	
Target 1	IL17RA
Target 1 Amino Acid Range	1 - 866
Target 1 Accession Number	NM_014339
Target 1 Description	Interleukin-17 receptor A
Target 1 Tag	EA
Target 2	IL17RC
Target 2 Amino Acid Range	1 - 720
Target 2 Accession Number	NM_153460
Target 2 Description	Interleukin-17 receptor C
Target 2 Tag	PK1
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete™ Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	IL-17 (IL-17A) (DiscoverX, 92-1229)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	10,000
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 (%)	98
Recovery After 48 Hours (%)	142
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 12.3
	Vial 2 13.9
	Vial 3 12.8
EC ₅₀ (g/mL)	Vial 1 16.2 x 10 ⁻⁹
	Vial 2 17.6 x 10 ⁻⁹
	Vial 3 16.9 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	March 2023
Expiration Date	March 2027

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: _____