

Certificate Of Analysis

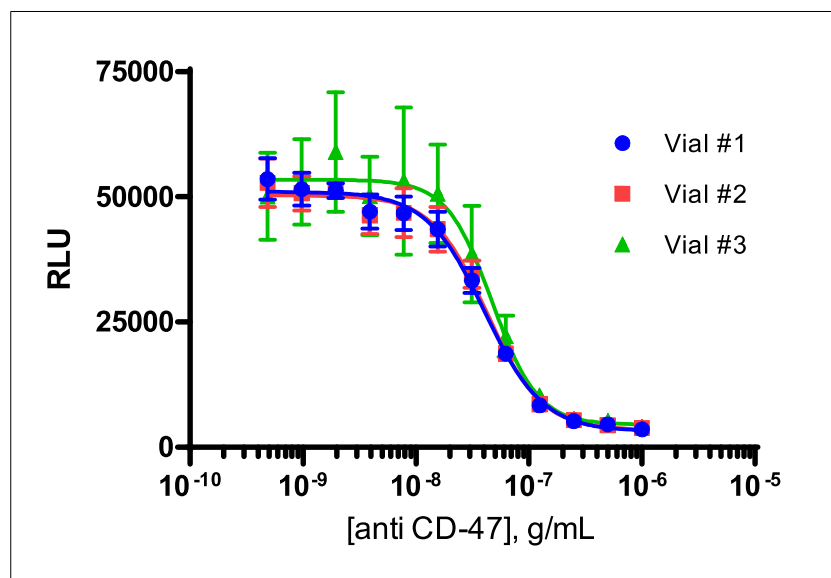
Background

Checkpoint Signaling assays are engineered to co-express a ProLink™ (PK) tagged immune checkpoint receptor and an Enzyme Acceptor (EA) tagged SH2 domain. Ligand engagement leads to receptor activation and phosphorylation, resulting in SH2-EA recruitment to the receptor, and forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

Product Name	PathHunter® Jurkat SIRPα Bioassay Cells
Cryovial Label	Jurkat SIRP α Checkpoint Signaling Bioassay Cells
Bioassay Catalog #	93-1135Y19
Bioassay Manufactured Lot #	22D2711
Passage # @ Freezing	9

Assay Information	
Target	SIRP α
Target Accession Number	NM_001040022.1
Target Description	Signal regulatory protein alpha
SH2 Domain	SHP1
Target Species	Human
Cell	Jurkat
CP Reagent	AssayComplete™ Cell Plating 0 Reagent (DiscoverX, 93-0563R0A)
Ligand	Jurkat CD47 Presenting Bioassay cells
Ligand Diluent	AssayComplete™ Cell Plating 0 Reagent
Detection Kit	PathHunter® Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	10,000
Cell Seeding Time (hours)	0
Ligand Incub Time (minutes)	1440
Agonist Incub Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	3.0
Fill Volume per Vial (mL)	0.2
Cell Viability	
Viability at Initial Thaw (%)	94
Recovery After 24 Hours (%)	134
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 15.0
	Vial 2 14.0
	Vial 3 12.5
IC ₅₀ (g/mL)	Vial 1 40.8 x 10 ⁻⁹
	Vial 2 42.7 x 10 ⁻⁹
	Vial 3 47.6 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	May 2022
Expiration Date	May 2025

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: _____