

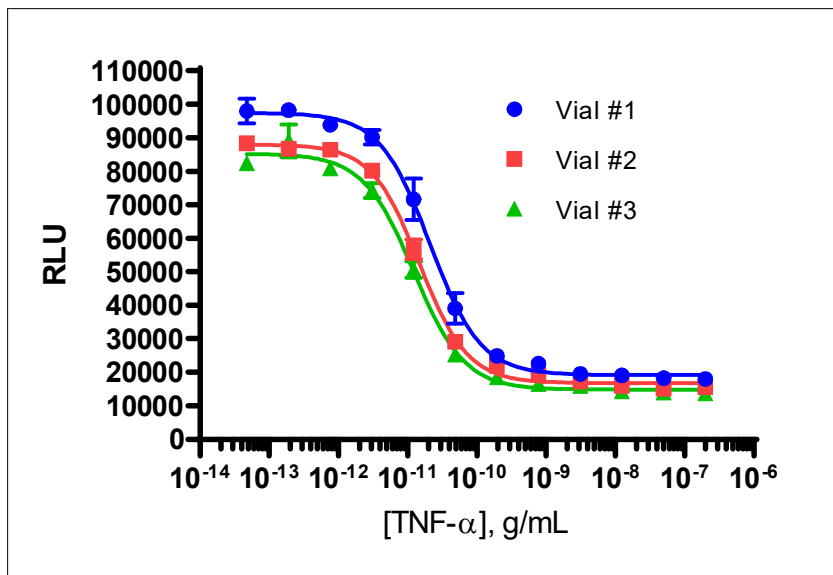
## Certificate Of Analysis

### Background

PathHunter® Degradation Bioassay cells are engineered to express an Enzyme Donor (ED) tagged target protein. Activation of the pathway results in degradation of the ED-tagged protein. Addition of exogenous Enzyme Acceptor (EA), and buffer, lyses the cell and forces complementation of the ED and EA enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes substrate to generate a chemiluminescent signal.

<b>Product Name</b>	<b>PathHunter® A549 IκB Bioassay Cells</b>
Cryovial Label	A549 IκB Degradation Bioassay Cells
Bioassay Catalog Number	93-0538B15
Bioassay Manufactured Lot #	22H1505
Passage # @ Freezing	4
<b>Assay Information</b>	
Target 1	IκB
Target 1 Accession Number	NM_020529
Target 1 Description	nuclear factor of kappa light polypeptide gene enhancer in B-cells inhibitor, alpha
Target 1 Tag	PL
Target 2	N/A
Target 2 Accession Number	N/A
Target 2 Description	N/A
Target 2 Tag	N/A
Target Species	Human
Cell	A549
CP Reagent	AssayComplete™ Cell Plating 38 Reagent (DiscoverX, 93-0563R38A)
Ligand	TNFα (human) (DiscoverX, 92-1097)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter® ED Bioassay Detection Kit (DiscoverX, 93-1043)
Cell Number/Well	10,000
Cell Seeding Time (hours)	48
Ligand Inc Time (minutes)	120
Agonist Inc Temperature (°C)	RT

Cell Density Information	
Cell Number (millions)	1.2
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	89
Recovery After 48 Hours (%)	277
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 5.4
	Vial 2 5.6
	Vial 3 6.0
EC <sub>50</sub> (g/mL)	Vial 1 0.021 x 10 <sup>-9</sup>
	Vial 2 0.015 x 10 <sup>-9</sup>
	Vial 3 0.012 x 10 <sup>-9</sup>



<b>Shipping and Storage Information</b>	
Shipping Conditions	Dry Ice
Storage Conditions	Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen.
Manufacturing Date	August 2022
Expiration Date	August 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: \_\_\_\_\_