

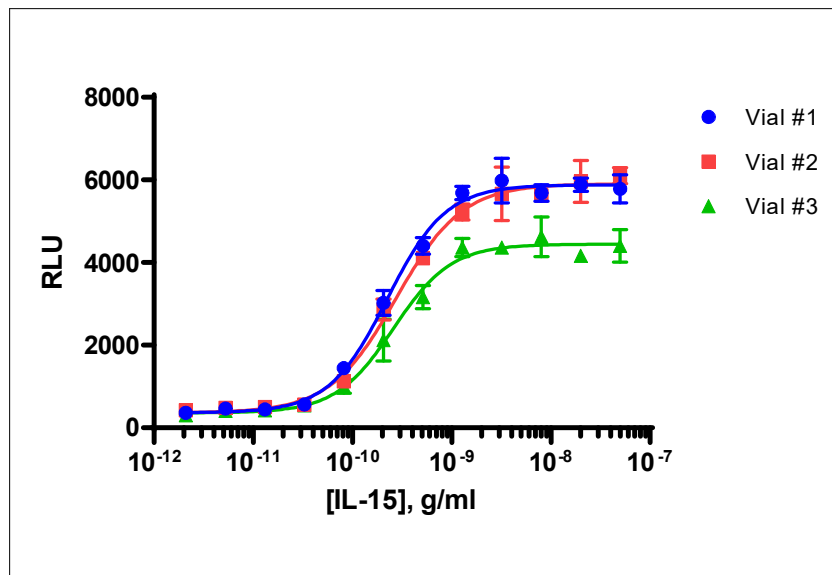
## Certificate Of Analysis

### Background

PathHunter Pathway Reporter Bioassay cells are engineered to express an Enzyme Donor (ED) tagged reporter protein controlled by a pathway-inducible transcriptional response element. Pathway activation via endogenous or exogenous target results in induced expressions 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® REH IL2RB-IL2RG STAT5 Signaling Pathway Reporter Cell Line</b>
Cryovial Label	IL-15 Reporter Bioassay Cells
Bioassay Catalog Number	93-1178Y055
Bioassay Manufactured Lot #	23K0208
Passage # @ Freezing	6
<b>Assay Information</b>	
Target 1	IL2RB
Target Species	Human
Target 1 Accession Number	NM_000878.3
Target 1 Description	Interleukin 2 receptor, beta
Reporter Response Element	STAT5
Reporter Tag	ePL
Cell	REH
CP Reagent	AssayComplete™ Cell Plating 0 Reagent (DiscoverX, 93-0563R0A)
Ligand	IL-15 (DiscoverX, 92-1265)
Ligand Diluent	AssayComplete™ Protein Dilution Buffer
Detection Kit	PathHunter® Detection Kit (DiscoverX, 93-0001)
Cell Number/Well	5000
Cell Seeding Time (hours)	0
Ligand Inc Time (minutes)	960
Agonist Inc Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	1.0
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	93
Recovery After 48 Hours (%)	137
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 16.1
	Vial 2 14.5
	Vial 3 14.7
EC <sub>50</sub> Check	Vial 1 2.2 x 10 <sup>-10</sup>
	Vial 2 2.7 x 10 <sup>-10</sup>
	Vial 3 2.6 x 10 <sup>-10</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	October 2023
Expiration Date	October 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: \_\_\_\_\_