

## Certificate Of Analysis

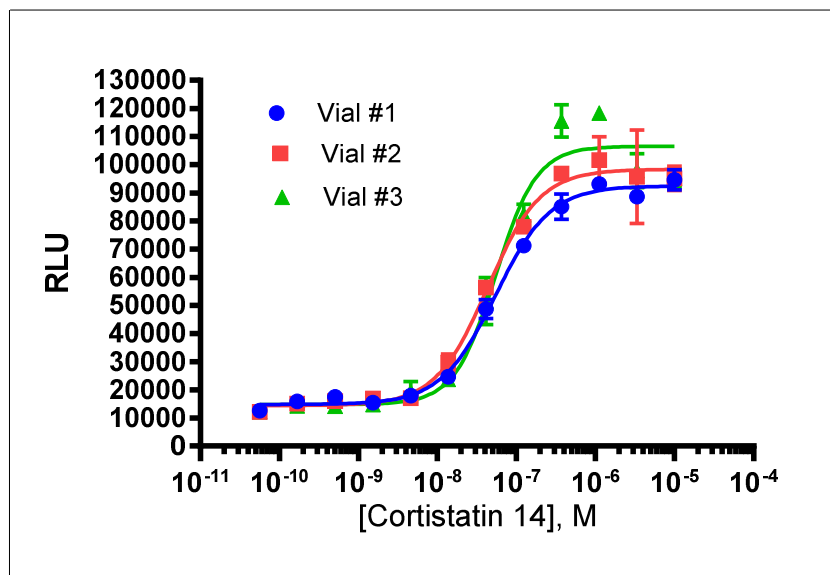
### Background

PathHunter<sup>®</sup>  $\beta$ -Arrestin GPCR Bioassay cells are engineered to co-express the ProLink<sup>™</sup> (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged  $\beta$ -Arrestin. Activation of the GPCR-PK induces  $\beta$ -Arrestin-EA recruitment, forcing complementation of the two  $\beta$ -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

<b>Product Name</b>	<b>PathHunter<sup>®</sup> CHO-K1 MRGPRX2 Bioassay Cells</b>
Cryovial Label	CHO-K1 MRGPRX2 Bioassay Cells
Bioassay Catalog #	93-0309Y2
Bioassay Manufactured Lot #	19D1901
Passage # @ Freezing	3

<b>Assay Information</b>	
Target 1	MRGPRX2
Target 1 Accession Number	NM_054030.2
Target 1 Description	MAS-related GPR, member X2
$\beta$ -Arrestin Isoform	$\beta$ -Arrestin-2
Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete <sup>™</sup> Cell Plating 2 Reagent (DiscoverX, 93-0563R2A)
Ligand	Cortistatin 14 (DiscoverX, 92-1076)
Ligand Diluent	AssayComplete <sup>™</sup> Cell Plating 2 Reagent
Detection Kit	PathHunter <sup>®</sup> Detection Kit (DiscoverX, 93-0001)
Cell Number/Well	5,000
Cell Seeding Time (hours)	24
Ligand Inc Time (minutes)	90
Agonist Inc Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	0.6
Fill Volume per Vial (mL)	0.1
Cell Viability	
Viability at Initial Thaw (%)	91
Recovery After 24 Hours (%)	114
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 7.5
	Vial 2 7.8
	Vial 3 7.7
EC <sub>50</sub> (M)	Vial 1 5.4 x 10 <sup>-8</sup>
	Vial 2 4.3 x 10 <sup>-8</sup>
	Vial 3 5.4 x 10 <sup>-8</sup>



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 2019
Expiration Date	April 2022

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: \_\_\_\_\_