

## Certificate Of Analysis

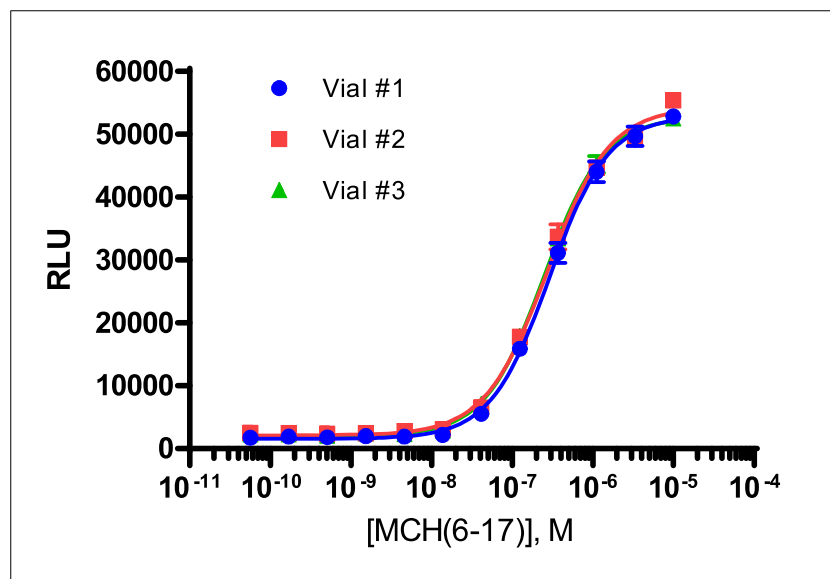
### Background

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> U2OS MCHR1 Bioassay Cells</b>
Cryovial Label	U2OS MCHR1 Beta-Arrestin Bioassay Cells
Bioassay Catalog #	93-0940Y3
Bioassay Manufactured Lot #	22J3001
Passage # @ Freezing	5

<b>Assay Information</b>	
Target 1	MCHR1
Target 1 Accession Number	AY562945
Target 1 Description	Melanin concentrating hormone receptor 1
$\beta$ -Arrestin Isoform	$\beta$ -Arrestin-2
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete <sup>™</sup> Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	Melanin Concentrating Hormone (6-17) (DiscoverX, 92-1173)
Ligand Diluent	AssayComplete <sup>™</sup> Cell Plating 5 Reagent
Detection Kit	PathHunter <sup>®</sup> Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	5,000
Cell Seeding Time (hours)	16
Ligand Inc Time (minutes)	90
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 (%)	92
Recovery After 48 Hours (%)	132
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 30.4
	Vial 2 22.0
	Vial 3 23.2
EC <sub>50</sub> (M)	Vial 1 288.4 x 10 <sup>-9</sup>
	Vial 2 271.3 x 10 <sup>-9</sup>
	Vial 3 245.8 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	October 2022
Expiration Date	October 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: \_\_\_\_\_