

Certificate Of Analysis

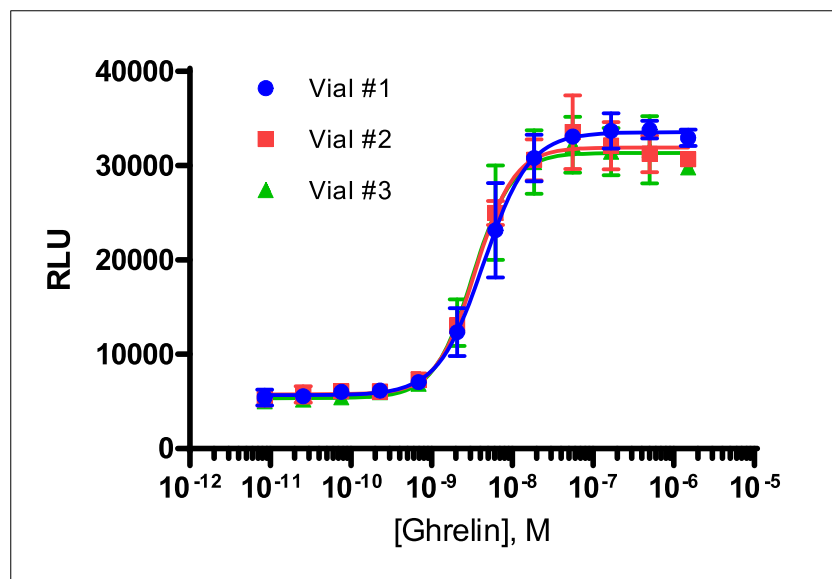
Background

PathHunter[®] β -Arrestin GPCR Bioassay cells are engineered to co-express the ProLink[™] (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged β -Arrestin. Activation of the GPCR-PK induces β -Arrestin-EA recruitment, 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[®] U2OS GHSR Bioassay Cells
Cryovial Label	U2OS GHSR Beta-Arrestin Bioassay Cells
Bioassay Catalog #	93-0242Y3
Bioassay Manufactured Lot #	22D1921
Passage # @ Freezing	3

Assay Information	
Target 1	GHSR
Target 1 Accession Number	NM_198407
Target 1 Description	Ghrelin receptor
β -Arrestin Isoform	β -Arrestin-2
Target Species	Human
Cell	U2OS
CP Reagent	AssayComplete [™] Cell Plating 5 Reagent (DiscoverX, 93-0563R5A)
Ligand	Ghrelin (DiscoverX, 92-1345)
Ligand Diluent	AssayComplete [™] Cell Plating 5 Reagent
Detection Kit	PathHunter [®] 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 (%)	90
Recovery After 48 Hours (%)	185
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 6.1
	Vial 2 5.9
	Vial 3 5.9
EC ₅₀ (M)	Vial 1 4.4 x 10 ⁻⁹
	Vial 2 3.5 x 10 ⁻⁹
	Vial 3 3.3 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	April 2022
Expiration Date	April 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: _____