

Certificate Of Analysis

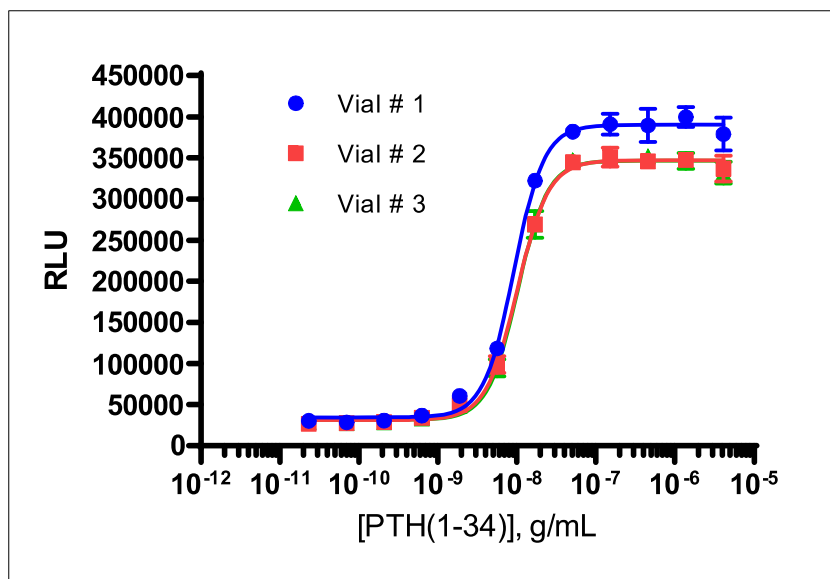
Background

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[®] CHO-K1 PTHR1 Bioassay
Cryovial Label	CHO-K1 PTHR1 Beta-Arrestin Bioassay Cells
Bioassay Catalog #	93-0315Y2
Bioassay Manufactured Lot #	22E3106
Passage # @ Freezing	3

Assay Information	
Target 1	PTHR1
Target 1 Accession Number	AY449732
Target 1 Description	Parathyroid hormone receptor 1
β -Arrestin Isoform	β -Arrestin-2
Target Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete [™] Cell Plating 0 Reagent (DiscoverX, 93-0563R0A)
Ligand	PTH(1-34) (DiscoverX, 92-1082)
Ligand Diluent	Protein Dilution Buffer
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933)
Cell Number/Well	10,000
Cell Seeding Time (hours)	48
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 (%)	96
Recovery After 24 Hours (%)	206
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 12.6
	Vial 2 12.8
	Vial 3 11.6
EC ₅₀ (g/mL)	Vial 1 9.2 x 10 ⁻⁹
	Vial 2 10.1 x 10 ⁻⁹
	Vial 3 10.3 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	June 2022
Expiration Date	June 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: _____