

CERTIFICATE OF ANALYSIS
Product Description and Information
BACKGROUND

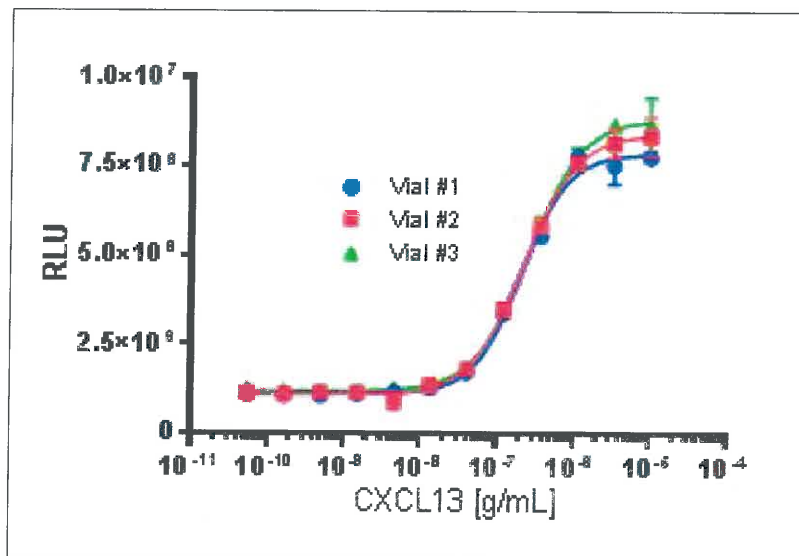
PathHunter[®] β -Arrestin GPCR Bioassay cells are engineered to co-express the ProLink[™] (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged β -Arrestin. Activation of the GPCR-PK induces β -Arrestin-EA recruitment, forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The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

Product Name	PathHunter [®] CHO-K1 CXCR5 Bioassay Cells
Cryovial Label	CHO-K1 CXCR5 Bioassay
Bulk Part #	93-0514Y2
Bulk Manufactured Lot #	15H2704
Bulk Passage # @ Freezing	3

Assay Information	
Target 1	CXCR5
Target 1 Accession Number	NM_001716
Target 1 Description	Burkitt lymphoma receptor
β -Arrestin Isoform	β -Arrestin-2
Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete [™] Cell Plating 2 Reagent (DiscoverRx, 93-0563R2A)
Ligand	CXCL13 (DiscoverRx, 92-1012)
Detection Kit	PathHunter [®] Bioassay Detection Kit (DiscoverRx, 93-0933)
Cell Number/Well	10000
Cell Seeding Time (hours)	24
Ligand Inc Time (minutes)	90
Agonist Inc Temperature (°C)	37

Product Specifications

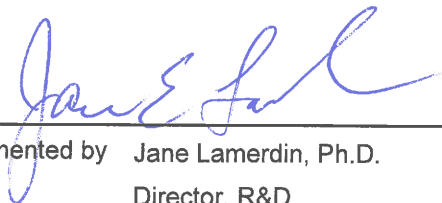
Cell Density Information		
Cell Number (millions)	1.25	
Fill Volume per Vial (mL)	0.1	
Cell Viability		
Viability at Initial Thaw (%)	96	
Recovery After 24 Hours (%)	95	
Mycoplasma and Sterility		
MycoSensor™ Mycoplasma Test	Passed	
Bacterial Sterility Test	Passed	
Functional Performance (3 manufactured vials)		
S:B Ratio	Vial 1	7.0
	Vial 2	7.4
	Vial 3	7.0
EC ₅₀ (g/mL)	Vial 1	2.09x10 ⁻⁷
	Vial 2	2.23x10 ⁻⁷
	Vial 3	2.54x10 ⁻⁷




Shipping and Storage Information

Shipping and Storage Information	
Shipping Conditions	Dry Ice
Storage Conditions	Short term (<2 weeks): -80°C; Long term (>2 weeks): Vapor phase of liquid N ₂ .
Date of Manufacture	August 2015
Expiration Date	August 2018

Shelf life of over 3 years has been established for PathHunter® cell lines in general when stored in the vapor phase of liquid nitrogen.

Signature: 
Documented by Jane Lamerdin, Ph.D.
Director, R&D

Date: 11/13/2015

Signature: 
Approved by Neil Charter, Ph.D.
Director, LeadHunter Services

Date: 11/13/2015