

## Certificate Of Analysis

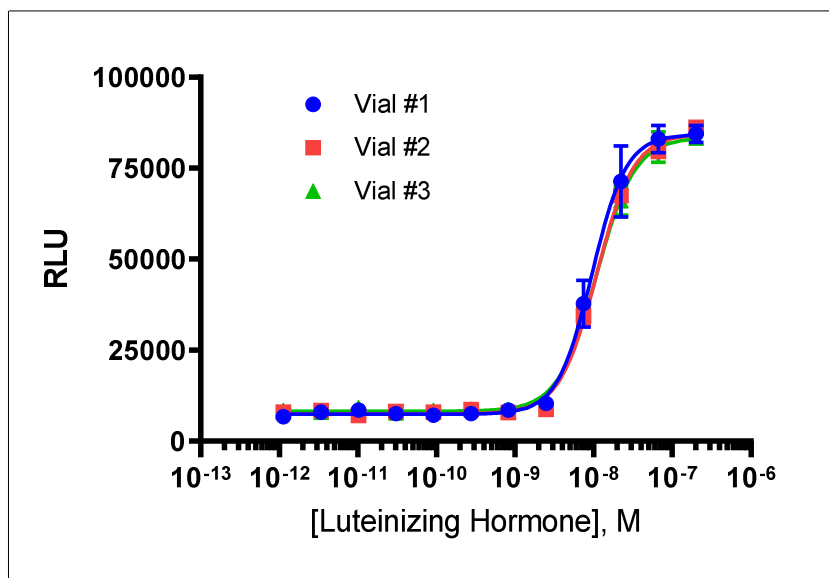
### Background

cAMP Hunter™ Bioassay cells are engineered to overexpress naturally Gi or Gs coupled, wildtype GPCRs and are designed to detect changes in intracellular cAMP levels in response to ligand stimulation of the receptor. These cells are designed to be used with the included HitHunter cAMP detection kit.

<b>Product Name</b>	<b>cAMP Hunter™ CHO-K1 LHCGR Bioassay Cells</b>
Cryovial Label	CHO-K1 LHCGR Bioassay Cells
Bioassay Catalog Number	95-0106Y2
Bioassay Manufactured Lot #	20B1005
Passage # @ Freezing	4

<b>Assay Information</b>	
Target	LHCGR
Accession Number	NM_000233
Target Description	Choriogonadotropin receptor
Receptor Family	Glycoprotein hormone receptor
Coupling	Gs
Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete™ Cell Plating 2 Reagent (DiscoverX, 93-0563R2A)
Ligand	Luteinizing hormone (DiscoverX, 92-1112)
Ligand Diluent	AssayComplete™ Protein Dilution Buffer B3
Detection Kit	HitHunter® cAMP Detection Kit (DiscoverX, 90-0075)
Cell Number/Well	30,000
Cell Seeding Time (hours)	48
Ligand Inc. Time (minutes)	30
Agonist Inc Temperature (°C)	37

Cell Density Information	
Cell Number (millions)	3.75
Fill Volume per Vial (mL)	0.2
Cell Viability	
Viability at Initial Thaw (%)	98
Recovery After 24 Hours (%)	136
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 12.6
	Vial 2 10.8
	Vial 3 9.8
EC <sub>50</sub> (M)	Vial 1 9.5 x 10 <sup>-9</sup>
	Vial 2 10.9 x 10 <sup>-9</sup>
	Vial 3 10.9 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	February 2020
Expiration Date	February 2023

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: \_\_\_\_\_