

Certificate Of Analysis

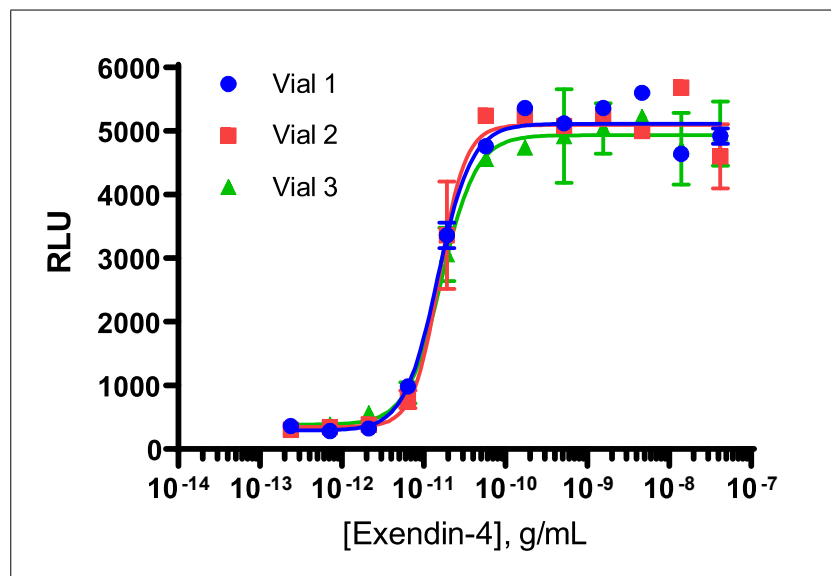
Background

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.

Product Name	cAMP Hunter™ CHO-K1 GLP1R Bioassay Cells
Cryovial Label	CHO-K1 GLP1R cAMP Bioassay Cells
Bioassay Catalog Number	95-0062Y2
Bioassay Manufactured Lot #	23G1911
Passage # @ Freezing	3

Assay Information	
Target	GLP1R
Accession Number	NM_002062.3
Target Description	Glucagon-like peptide receptor 1
Receptor Family	Glucagon
Coupling	Gs
Target Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete™ Cell Plating 2 Reagent (DiscoverX, 93-0563R2A)
Ligand	Exendin-4 (DiscoverX, 92-1115)
Ligand Diluent	AssayComplete™ Protein Dilution Buffer B2
Detection Kit	HitHunter® cAMP Detection Kit (DiscoverX, 90-0075)
Cell Number/Well	30,000
Cell Seeding Time (hours)	24
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 (%)	213
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 13.7
	Vial 2 15.3
	Vial 3 14.6
EC ₅₀ (g/mL)	Vial 1 0.015 x 10 ⁻⁹
	Vial 2 0.016 x 10 ⁻⁹
	Vial 3 0.016 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	July 2023
Expiration Date	July 2026

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: _____