

## Certificate Of Analysis

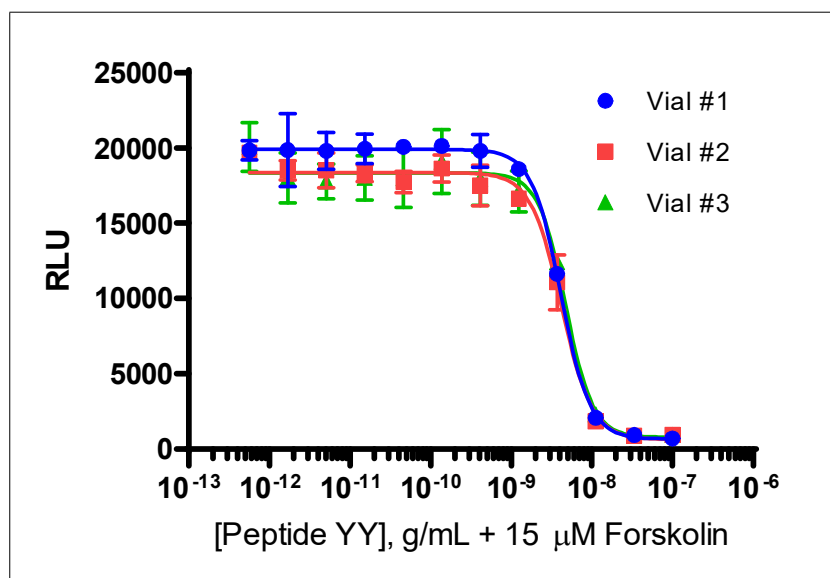
### Background

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 NPY2R Bioassay Cells</b>
Cryovial Label	CHO-K1 NPY2R cAMP Bioassay Cells
Bioassay Catalog Number	95-0077Y2
Bioassay Manufactured Lot #	22B2808
Passage # @ Freezing	3

<b>Assay Information</b>	
Target	NPY2R
Accession Number	AY236540
Target Description	Neuropeptide Y receptor Y2
Receptor Family	Neuropeptide Y
Coupling	Gi/Go & Gq
Target Species	Human
Cell	CHO-K1
CP Reagent	AssayComplete™ Cell Plating 2 Reagent (DiscoverX, 93-0563R2A)
Ligand	Peptide YY, Human (DiscoverX, 92-1063)
Ligand Diluent	AssayComplete™ cAMP Hunter Cell Assay Buffer
Detection Kit	cAMP Detection Kit for Bioassays (DiscoverX, 90-0075B2)
Cell Number/Well	10,000
Cell Seeding Time (hours)	48
Ligand Inc. Time (minutes)	30
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 (%)	94
Recovery After 24 Hours (%)	157
Mycoplasma and Sterility	
Mycoplasma Test	Passed
Sterility Test	Passed
Functional Performance (3 manufactured vials)	
S:B Ratio	Vial 1 28.3
	Vial 2 21.4
	Vial 3 21.4
EC <sub>50</sub> (g/mL)	Vial 1 4.1 x 10 <sup>-9</sup>
	Vial 2 4.2 x 10 <sup>-9</sup>
	Vial 3 4.7 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	March 2022
Expiration Date	March 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: \_\_\_\_\_