

cAMP Hunter™ U2OS mGLP1R Gs Cell Line**Catalog Number:** 95-0179C3**Lot Number:**

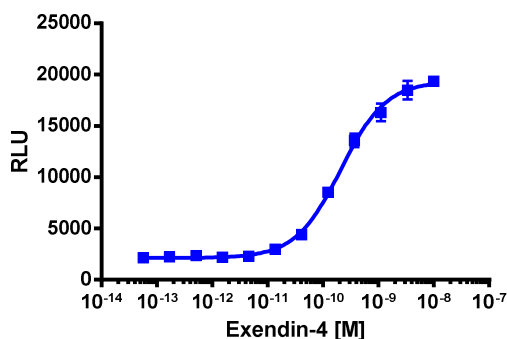
See Vial

Contents: 2 vials, 1 x 10⁶ cells per vial in 1 mL**Background**

cAMP Hunter™ Gs cell lines overexpress naturally Gs coupled, wildtype GPCRs and are designed to detect increases in intracellular cAMP levels in response to agonist stimulation of the receptor. These cell lines are designed to be used in conjunction with the HitHunter® cAMP Assay Detection Kit.

Product Information**Target GPCR:** mGLP1R**Description:** Glucagon-like peptide receptor 1**Receptor Family:** Glucagon**Coupling:** Gs**Accession Number:** NM_021332**GPCR Species:** Mouse**Cell Type:** U2OS**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.**Functional Performance**

Cells were plated in a 384-well plate and incubated overnight at 37°C and 5% CO₂ to allow the cells to attach and grow. Cells were then stimulated with a control agonist, using the assay conditions described below. Following stimulation, signal was detected using the HitHunter cAMP Assay Detection Kit according to the recommended protocol. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.

**Cell Number/Well:**

10000

Control Agonist:

Exendin-4

Agonist Incubation Time (minutes):

30

Agonist Incubation Temperature (°C):

37

EC₅₀ for Agonist Stimulation (nM):

0.2

Signal:Background at Agonist E_{max}:

9.1

Important! This assay requires an additional step: Please refer to Additional Protocol Information section.

Passage Stability

This cell line has been confirmed to be stable through 10 passages with no significant drop in assay window or change in EC₅₀.

Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	HitHunter® cAMP Assay for Small Molecules HitHunter cAMP Assay for Biologics	90-0075SM Series 90-0075LM Series
Cell Culture	AssayComplete™ Cell Culture Kit-103	92-3103G
Cell Plating	AssayComplete™ Cell Plating 5 Reagent	93-0563R5A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T3	92-4103TR
Cell Freezing	AssayComplete™ Freezing Reagent F3	92-5103FR

*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	0.25	92-0028
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	500	92-0030

Additional Ligand Information

Control Agonist: Exendin-4

Vendor: DiscoverX® (Catalog No. 92-1115)

Additional Protocol Information

Additional Cell Line Info: Cell line also over-expresses ADCY2 (NM_020546.2) to enhance stimulation of cAMP synthesis in response to activation of the Gs-coupled GPCR.

Additional Protocol Notes: It is recommended that the control compound be serially diluted in a buffer (HBSS/10 mM HEPES or PBS) + 1% FBS to ensure optimal assay performance.

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail info@discoverx.com. For additional information, please visit discoverx.com.

Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit discoverx.com/license for a list of products that are governed by limited use label license terms and relevant patent and trademark information.