

## cAMP Hunter™ CHO-K1 GIPR Gs Cell Line

**Catalog Number:** 95-0146C2

**Lot Number:** See Vial

**Contents:** 2 x 10<sup>6</sup> cells per vial in 1 mL

### Background

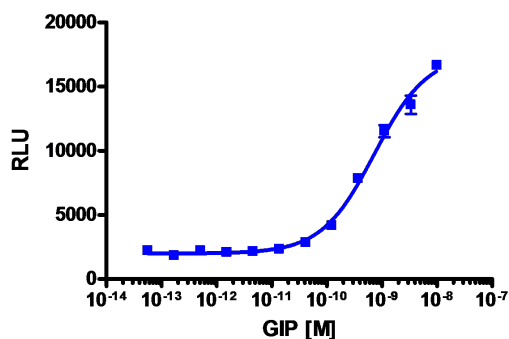
cAMP Hunter™ Gs cell lines overexpress naturally Gs coupled, wildtype GPCRs and are designed to detect increases in intracellular cAMP levels in response to agonist stimulation of the receptor. These cell lines are designed to be used in conjunction with the HitHunter® cAMP Assay Detection Kit.

### Product Information

**Target GPCR:** GIPR  
**Description:** Gastric inhibitory polypeptide receptor  
**Receptor Family:** Glucagon  
**Coupling:** Gs  
**Accession Number:** NM\_000164  
**GPCR Species:** Human  
**Cell Type:** CHO-K1  
**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% CO<sub>2</sub> 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:** GIP (Gastric Inhibitory Peptide), Human  
**Agonist Incubation Time (minutes):** 30  
**Agonist Incubation Temperature (°C):** 37  
**EC<sub>50</sub> for Agonist Stimulation (nM):** 0.73  
**Signal:Background at Agonist E<sub>max</sub>:** 7.5

### Passage Stability

This cell line has been confirmed to be stable through a minimum of 10 passages with no significant drop in assay window or change in EC<sub>50</sub>.

### 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 <sup>®</sup> cAMP Assay for Small Molecules HitHunter cAMP Assay for Biologics	90-0075SM Series 90-0075LM Series
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-107	92-3107G
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 2 Reagent	93-0563R2A
Cell Detachment	AssayComplete <sup>™</sup> Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T2	92-4102TR
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F2	92-5102FR

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

### Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete <sup>™</sup> Puromycin	Not Applicable	Not Applicable
AssayComplete <sup>™</sup> Hygromycin B	Not Applicable	Not Applicable
AssayComplete <sup>™</sup> G418	800	92-0030

### Additional Ligand Information

**Control Agonist:** GIP (Gastric Inhibitory Peptide), Human

**Vendor:** Eurofins DiscoverX<sup>®</sup> (Catalog No. 92-1078)

### 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 <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.