

cAMP Hunter™ CHO-K1 GPR35 Gi Cell Line

Catalog Number: 95-0152C2 **Lot Number:** See Vial
Contents: 2 vials, 1 x 10⁶ cells per vial in 1 mL

Background

cAMP Hunter™ Gi cell lines overexpress naturally Gi coupled, wildtype GPCRs and are designed to detect inhibition of intracellular cAMP production 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: GPR35
Description: G-protein coupled receptor 35
Receptor Family: Class A Orphan
Coupling: Gi/Go
Accession Number: NM_005301
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% CO₂ to allow the cells to attach and grow. Cells were then stimulated with a control agonist + forskolin (Figure 1A) or with forskolin only (Figure 1B), 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: 5000
Control Agonist: Zaprinast
Agonist Incubation Time (minutes): 30
Agonist Incubation Temperature (°C): 37

Figure 1A. Agonist Dose Response

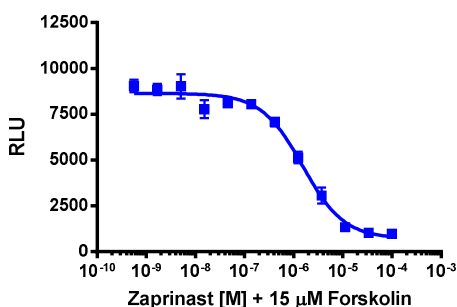
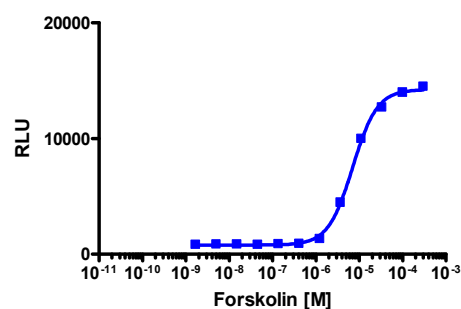


Figure 1B. Forskolin Dose Response



EC₅₀ for Agonist Stimulation (nM): 1520.5
Signal:Background at Agonist E_{max}: 9.2
Forskolin Concentration (μM): 15

EC₅₀ for Forskolin Stimulation (μM): 7.0
Signal:Background at Forskolin E_{max}: 18.0

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

*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	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	800	92-0030

Additional Ligand Information

Control Agonist: Zaprinast

Vendor: DiscoverX® (Catalog No. 92-1048)

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.