

## cAMP Hunter™ CHO-K1 HTR7A Gs Cell Line

**Catalog Number:** 95-0151C2 **Lot Number:** See Vial  
**Contents:** 2 vials, 1 x 10<sup>6</sup> cells per vial in 1 mL

### Background

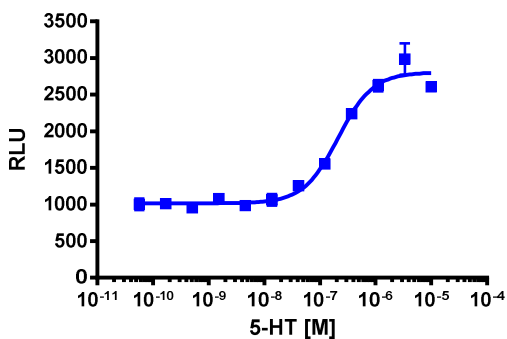
cAMP Hunter™ Gs cell lines overexpress naturally Gs coupled, wildtype GPCRs and are designed to detect increases in intracellular cAMP levels in response to agonist stimulation of the receptor. These cell lines are designed to be used in conjunction with the HitHunter® cAMP Assay Detection Kit.

### Product Information

**Target GPCR:** HTR7A  
**Description:** 5-hydroxytryptamine receptor 7A  
**Receptor Family:** 5-Hydroxytryptamine  
**Coupling:** Gs  
**Accession Number:** NM\_000872.4  
**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:** 5-HT  
**Agonist Incubation Time (minutes):** 30  
**Agonist Incubation Temperature (°C):** 37  
**EC<sub>50</sub> for Agonist Stimulation (nM):** 212  
**Signal:Background at Agonist E<sub>max</sub>:** 3

### Passage Stability

This cell line has been confirmed to be stable through 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® cAMP Assay for Small Molecules<br>HitHunter cAMP Assay for Biologics | 90-0075SM Series<br>90-0075LM Series |
| Cell Culture    | AssayComplete™ Cell Culture Kit-108   | 92-3108G                             |
| Cell Plating    | AssayComplete™ Cell Plating 21 Reagent  | 93-0563R21A                          |
| 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:** 5-HT

**Vendor:** DiscoverX® (Catalog No. 92-1108)

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](mailto:info@discoverx.com). For additional information, please visit [discoverx.com](http://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](https://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.