



GIP Receptor (human) Reporter Assay Kit

Item No. 502875

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

| | | |
|-----------------------------|----|---|
| GENERAL INFORMATION | 3 | Materials Supplied |
| | 3 | Safety Data |
| | 4 | Precautions |
| | 4 | If You Have Problems |
| | 5 | Storage and Stability |
| | 5 | Materials Needed but Not Supplied |
| INTRODUCTION | 6 | Background |
| | 7 | About This Assay |
| | 8 | Principle Of This Assay |
| ASSAY PROTOCOL | 9 | Plate Set Up |
| | 10 | Addition of Cells to the Reverse Transfection Plate |
| | 11 | Cell Stimulation |
| | 12 | Performing the SEAP Assay |
| PERFORMANCE CHARACTERISTICS | 13 | Calculations |
| | 14 | Performance Characteristics |
| RESOURCES | 15 | Troubleshooting |
| | 16 | References |
| | 17 | Plate Template |
| | 18 | Notes |
| | 19 | Warranty and Limitation of Remedy |

GENERAL INFORMATION

Materials Supplied

This kit will arrive packaged as a -20°C kit. After opening the kit, store individual components as stated below.

| Item Number | Item | Quantity/ Size | Storage |
|-------------|---|-------------------|---------|
| 401126 | GIP Receptor (human) Reverse Transfection Strip Plate | 1 plate | -20°C |
| 401127 | Tirzepatide Positive Control (0.1 mM) | 1 vial/25 µl | -20°C |
| 600183 | SEAP Substrate (Luminescence) | 1 vial/15 ml | 4°C |
| 700029 | 96-Well Solid Plate (white) | 3 plates | RT |
| 400012 | 96-Well Cover Sheet | 3 ea | RT |

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



WARNING: THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

Safety Data

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent *via* email to your institution.

Precautions

Please read these instructions carefully before beginning this assay.

The reagents in this kit have been tested and formulated to work exclusively with Cayman Chemical's GIP Receptor (human) Reporter Assay Kit. This kit may not perform as described if any reagent or procedure is replaced or modified.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888

Email: techserv@caymanchem.com

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored as directed in the Materials Supplied section (see page 3) and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. HEK293T (or HEK293T/17) cells, available from ATCC
2. Dulbecco's Modified Eagle's Medium (DMEM)
3. Fetal bovine serum (FBS)
4. Penicillin-streptomycin (100X) (ThermoFisher™ Scientific, Catalog No. 15140-122)
5. A plate reader capable of measuring luminescence
6. Adjustable and multichannel pipettes with pipette tips
7. An incubator/oven set at 65°C

Background

Glucose-dependent insulintropic polypeptide receptor (GIPR) is a transmembrane glycoprotein hormone receptor and class B1 G protein-coupled receptor (GPCR).¹ It is expressed in adipose tissue and the pancreas, gut, heart, pituitary gland, adrenal cortex, and brain.^{2,3} GIPR is activated by the endogenous ligands GIP (1-42) and GIP (1-30) amide, inducing primarily $G\alpha_s$ signaling, to regulate insulin and glucagon secretion, bone remodeling, and lipid metabolism.^{2,4,5} Knockout of *GIPR* increases lipolysis in white adipose tissue, diminishes body weight gain and reduces fat mass in a mouse model of obesity induced by a high-fat diet, but both pharmacologic agonism and antagonism of GIPR induces weight loss and improves metabolic outcomes in preclinical models of obesity and diabetes.^{4,6} Pharmacologic agonism of GIPR is associated with a decreased risk of the development of non-alcoholic fatty liver disease (NAFLD).⁷ These effects highlight the importance of continued research and development of GIPR agonists as potential therapeutics for diabetes and obesity.

About This Assay

Cayman's reverse transfection reporter assays have overcome many of the disadvantages of other transfection approaches. In this method, a proprietary transfection complex containing DNA and an optimized mixture of lipids and proteins has been evenly applied and processed on the culture surface of multi-well plates. Adherent cells, supplied by the user, are applied directly to the plate and allowed to grow in the coated wells. Using this method, the uptake of the DNA complex by the cell increases dramatically compared to solution-phase transfection, enhancing both the transfection efficiency and co-transfection efficiency for multiple plasmids.

Cayman's GIP Receptor (human) Reporter Assay Kit consists of a 96-well plate coated with a transfection complex containing DNA constructs for expressing GIPR and a cAMP response element-regulated secreted alkaline phosphatase (SEAP) reporter (GIP Receptor Reverse Transfection Strip Plate). Cells grown on the transfection complex will express GIPR at the cell surface within 24 hours. Binding of agonists to GIPR initiates a signal transduction cascade through the $G\alpha_s$ -coupled adenylate cyclase pathway resulting in the expression of SEAP, which is secreted into the cell culture medium. Aliquots of culture medium are collected at 6-8 hours after stimulation and SEAP activity is measured following the addition of a luminescence-based alkaline phosphatase substrate provided in the kit (SEAP Substrate). The kit is easy to use and can be readily applied to high-throughput screening for therapeutic compounds regulating the activation of GIPR. A synthetic peptide agonist, tirzepatide, is included for use as a positive control. The kit provides a sufficient volume of reagents to measure SEAP activity at three time points using the three included white assay plates.

Principle Of This Assay

The binding of an agonist to GIPR activates $G\alpha_s$ and triggers the cAMP/PKA/ CREB signaling pathway. Phosphorylated CREB binds to the CREB-responsive element (CRE) in the promoter region of the co-transfected reporter, which induces the expression of SEAP. The activity of SEAP is measured as a luminescence signal produced after incubation with the provided substrate.

ASSAY PROTOCOL

Plate Set Up

There is no specific pattern for using the wells on the plate. A typical experiment will include wells containing cells treated with tirzepatide as a positive control, cells treated with experimental compounds, and untreated cells. It is recommended that each treatment be performed at least in triplicate. To determine the EC_{50} value of a compound, several concentrations of the compound should be included in the assay. The volume of Tirzepatide Positive Control (0.1 mM) is sufficient to run a full dose-response curve with replicates or to serve as a control at approximately EC_{80} for testing inhibitors on the whole plate. Record the contents of each well on the template sheet provided on page 17.

Addition of Cells to the Reverse Transfection Plate

IMPORTANT

Before starting the experiment, pre-warm the required volume of culture medium and make sure a sufficient number of actively growing cells are available.

1. Remove the GIP Receptor (human) Reverse Transfection Strip Plate (Item No. 401126) from the freezer and allow it to equilibrate to room temperature inside the sealed bag. After it has reached room temperature, clean the bag with 70% alcohol before opening it and place the plate inside a cell culture hood.

NOTE: If not using the whole plate within one experiment, remove the number of strips needed, place the remaining strips back in the bag, and store them in a desiccator protected from UV light at room temperature for up to a week. Alternatively, the remaining strips can be sealed in the bag with the desiccant pack and stored at -20°C for up to two months.

2. Seed HEK293T cells at a density of 40,000-60,000 cells/well in 200 µl of DMEM containing 10% FBS and 1X penicillin/streptomycin. Leave the plate in the hood for 30-45 minutes.
3. Place the plate in a 37°C CO₂ incubator and incubate for 16-20 hours.

Cell Stimulation

1. After 16-20 hours of incubation, carefully aspirate the culture medium from each well.
2. Replenish the cells with 150 µl of pre-warmed stimulation medium (DMEM with 0.5% FBS) per well.
3. Prepare test compounds at 4X the desired final concentration in DMEM and pipette 50 µL into the assigned wells. Pipette 50 µL of DMEM into wells designated for untreated cells. For positive control wells, dilute the Tirzepatide Positive Control (0.1 mM) (Item No. 401127) 1:250 in serum-free DMEM and add 50 µl to the positive control wells. At this concentration (100 nM), tirzepatide will induce a >5-fold increase in SEAP activity in 6-8 hours compared to the untreated control.

Pipetting Hints

- Use different tips to pipette each reagent.
- Avoid introducing bubbles to the well.
- Do not expose the pipette tip to the reagent(s) already in the well.

Before performing the assay, remove the SEAP Substrate (Luminescence) (Item No. 600183) from the refrigerator and allow it to equilibrate to room temperature.

1. After 6-8 hours of stimulation with test compounds and controls, transfer the plate from the incubator to a cell culture hood.
2. Inside the culture hood, use a multi-channel pipette to gently pipette up and down a few times. Collect 10 µl of medium from each well and transfer it to the corresponding well of a 96-Well Solid Plate (white) (Item No. 700029).

NOTE: Avoid contact of the pipette tip with the plate bottom to minimize disruption of the cell layer. Perform this step inside the cell culture hood and return the plate to the incubator if sampling at later time point(s) is needed.

3. Cover the assay plate with the 96-Well Cover Sheet (Item No. 400012) and incubate the plate in a 65°C incubator for 30 minutes to heat inactivate endogenous alkaline phosphatase. The SEAP expressed in this assay is stable under this condition.
4. Remove the plate from the 65°C incubator, remove the cover sheet, and allow the plate to cool to room temperature.
5. Add 50 µl SEAP Substrate to each well, shake/tap briefly to mix, and incubate the plate at room temperature for 5-15 minutes.
6. Scan the plate for luminescence in a microplate reader.

NOTE: The plate should be read immediately after 5-15 minutes of incubation with SEAP Substrate. When multiple plates are processed at the same time, the time interval between plates for the addition of substrate and plate reading should be consistent.

Calculations

Determination of EC₅₀

The term half-maximal effective concentration (EC₅₀) refers to the concentration of a drug that induces a response halfway between the baseline and maximum after a specific exposure time. The dose-response curve of a typical agonist follows a sigmoidal curve with a bottom plateau (untreated cells) and a top plateau (drug saturation). See Figure 1 on page 14 for a typical GIPR agonist dose-response curve.

For each compound, normalize the relative luminescent unit (RLU) results to run from 0% (no drug added) to 100% (saturating dose) by using the following formula:

% Response at X Concentration =

$$\left[\frac{(\text{RLU at X concentration}) - (\text{RLU of untreated cells})}{\text{Maximal RLU (saturation)} - (\text{RLU of untreated cells})} \right] \times 100$$

Graph the percent response versus log drug concentration. In the resulting sigmoidal dose-response curve, find the best-fit value for the log EC₅₀ (the concentration that gives a 50% response; the middle of the curve).

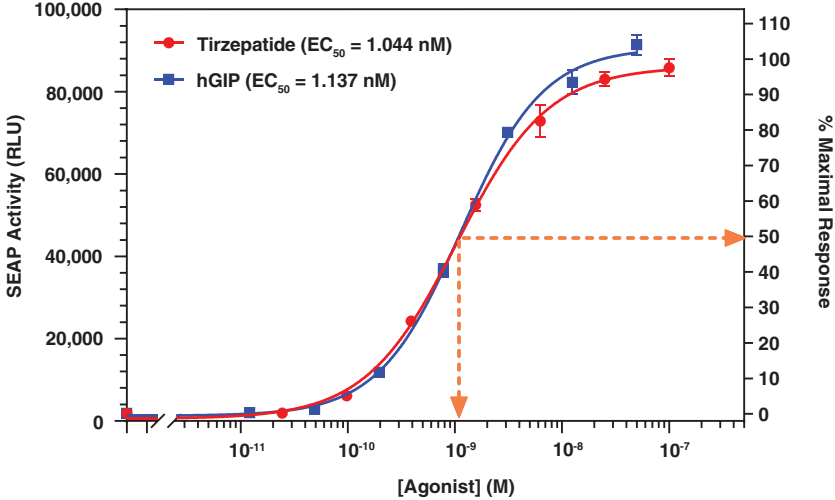


Figure 1. SEAP activity in HEK293T cells transiently-transfected with GIPR in response to human GIP and tirzepatide stimulation. HEK293T cells were seeded on a GIP receptor (human) Reverse Transfection Strip Plate at a density of 40,000 cells/well and incubated overnight. The next day, the cells were treated with different doses of human GIP (hGIP) up to 50 nM or tirzepatide up to 100 nM in reduced serum medium. After six hours of stimulation, 10 µl of culture medium was collected from each well and the SEAP activity of each sample was measured according to the Performing the SEAP Assay section (see page 12). The calculated EC₅₀ values for hGIP and tirzepatide from the fitted curves are 1.137 nM and 1.044 nM, respectively. The Z' value of the assay is >0.9.

NOTE: The fold of stimulation, Z' value, and calculated EC₅₀ may vary with cell lines, cell passage, and culture conditions.

Troubleshooting

| Problem | Possible Causes | Recommended Solutions |
|--|--|--|
| Dispersion of replicates or erratic response curve of test compounds | A. Uneven cell distribution B. Poor pipetting C. Not well mixed when sampling D. Bubble in assay wells | A. Make sure cells are in homogenous suspension at plating and allow the cells to sit for 30-45 min before putting into incubator B. Pipette carefully C. Pipette up and down a few times before collecting sample D. Carefully tap the side of the plate to remove bubbles |
| Low reading in wells | A. Reading time is too short B. Samples overheated/dried C. The substrate is too cold | A. Increase the integration time B. Do not heat-inactivate the sample plate above 65°C or longer than 30 minutes C. Warm up the SEAP substrate to room temperature before use |
| Sample signal is too strong | A. Cell density was too high B. Insufficient heat inactivation of endogenous alkaline phosphatase activity | A. Reduce cell plating density B. Correct the duration or temperature of heat inactivation step |
| Poor control curve/signal | A. Control compound degraded B. Pipetting error C. Splashing of sample D. Volume carry-over during dilution | A. Avoid free-thaw of positive control B. Check pipette volume C. Dispense carefully D. Use new tip for each pipetting |

References

1. Aksu, H., Demirbilek, A., and Uba, A.I. Insights into the structure and activation mechanism of some class B1 GPCR family members. *Mol. Biol. Rep.* **51**(1), 966 (2024).

2. Nauck, M.A., Quast, D.R., Wefers, J., *et al.* The evolving story of incretins (GIP and GLP-1) in metabolic and cardiovascular disease: A pathophysiological update. *Diabetes Obes. Metab.* **23**(Suppl 3), 5-29 (2021).

3. Whitaker, G.M., Lynn, F.C., McIntosh, C.H.S., *et al.* Regulation of GIP and GLP1 receptor cell surface expression by N-glycosylation and receptor heteromerization. *PLoS One* **7**(3), e32675 (2012).

4. Boer, G.A., Keenan, S.N., Miotto, P.M., *et al.* GIP receptor deletion in mice confers resistance to high-fat diet-induced obesity via alterations in energy expenditure and adipose tissue lipid metabolism. *Am. J. Physiol. Endocrinol. Metab.* **320**(4), E835-E845 (2021).

5. Rosenkilde, M.M., Lindquist, P., Kizilkaya, H.S., *et al.* GIP-derived GIP receptor antagonists - a review of their role in GIP receptor pharmacology. *Peptides* **177**, 171212 (2024).

6. James-Okoro, P.-P., Lewis, E. C., Gribble, F.M., *et al.* The role of GIPR in food intake control. *Front. Endocrinol.* **16**, 1532076 (2025).

7. Xie, S., Chen, D., Cai, Y., *et al.* Evaluating the efficacy of GIPR agonists on non-alcoholic fatty liver disease: A mediation mendelian randomization study. *Digestive and Liver Disease* **56**(10), 1730-1737 (2024).

| | | | | | | | | |
|----|---|---|---|---|---|---|---|---|
| 12 | | | | | | | | |
| 11 | | | | | | | | |
| 10 | | | | | | | | |
| 9 | | | | | | | | |
| 8 | | | | | | | | |
| 7 | | | | | | | | |
| 6 | | | | | | | | |
| 5 | | | | | | | | |
| 4 | | | | | | | | |
| 3 | | | | | | | | |
| 2 | | | | | | | | |
| 1 | | | | | | | | |
| | A | B | C | D | E | F | G | H |

Warranty and Limitation of Remedy

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©08/14/2025, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.
Printed in U.S.A.

