

**Description**

Rat GLP-1R/CRE Luciferase Reporter HEK293 Cell Line is an engineered HEK293 cell line expressing firefly luciferase under the control of cAMP response element (CRE), and rat GLP-1R (*Rattus norvegicus* Glucagon-like peptide 1 receptor; NM\_012728.2). Activation of rat GLP-1R in these cells can be monitored by measuring luciferase activity.

The functionality of Rat GLP-1R/CRE Luciferase Reporter HEK293 Cell Line was validated in dose-response assays using the peptide agonists human Glucagon-like peptide 1 (7-37).

Rat GLP-1R/CRE Luciferase Reporter HEK293 Cell Line is a *Rattus norvegicus* version of GLP-1R/CRE Luciferase Reporter HEK293 Cell Line (BPS Bioscience #78176).

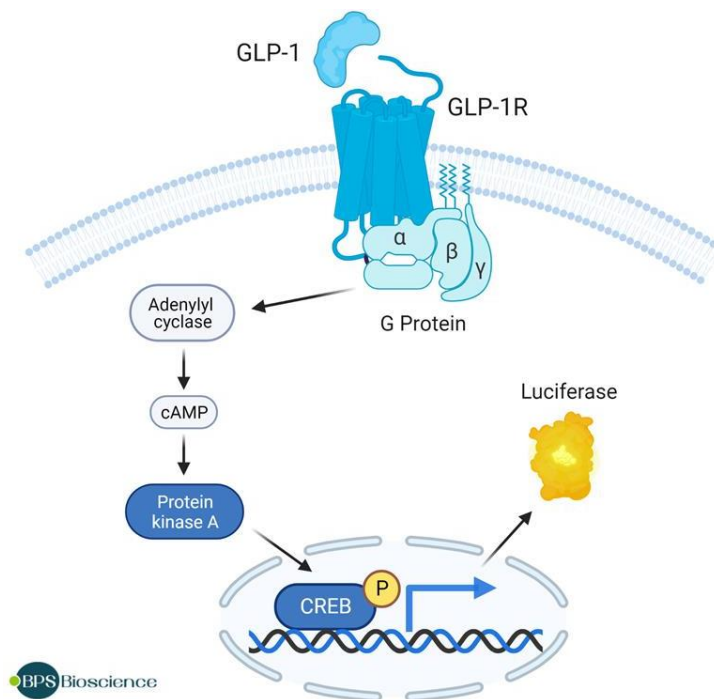


Figure 1. Illustration of mechanism of action in Rat GLP-1R/CRE Luciferase Reporter HEK293 Cell Line.

**Background**

GLP-1R, a member of the class B family of G protein-coupled receptors (GPCRs or secretin-like receptors), is a transmembrane protein primarily found in pancreatic  $\beta$  cells and brain neurons. GLP-1R is activated by the peptide hormone glucagon-like peptide 1 (GLP-1), which has two active forms, GLP-1 (7-37) and GLP-1 (7-36) amide. GLP-1R plays an important role in controlling blood sugar level by enhancing glucose-stimulated insulin secretion, glucose, lipid metabolism, and satiety. Its role in the brain seems to be in the control of appetite. Research efforts have focused on the regulation of the GLP-1R mediated signaling pathway as a therapeutic approach to type 2 diabetes (T2DM) and have resulted in the development of several GLP-1 FDA-approved agonists. In addition to their role in insulin secretion, GLP-1R agonists can also contribute to weight management, decrease the potential for cardiovascular diseases and protect beta cells. A role in tumor development in patients with T2DM is also being investigated but further studies are required to fully understand the functions of GLP-1R and its agonists.

**Application**

Screen for agonists of rat GLP-1R in a cellular model.

**Materials Provided**

Components	Format
2 vials of frozen cells	Each vial contains >1 x 10 <sup>6</sup> cells in 1 ml of Cell Freezing Medium (BPS Bioscience #79796)

**Host Cell**

HEK293, Human Embryonic Kidney, epithelial-like cells, adherent.

**Mycoplasma Testing**

The cell line has been screened to confirm the absence of Mycoplasma species.

**Materials Required but Not Supplied**

These materials are not supplied with this cell line but are necessary for cell culture and cellular assays. BPS Bioscience reagents systems are validated and optimized for use with this cell line and are highly recommended for best results. Media components are provided in the Media Formulations section.

*Media Required for Cell Culture*

Name	Ordering Information
Thaw Medium 1	BPS Bioscience #60187
Growth Medium 1U	BPS Bioscience #78548

*Materials Required for Cellular Assay*

Name	Ordering Information
GLP-1 (7-37)	BPS Bioscience #82667
Opti-MEM reduced serum medium (Assay Medium)	Thermo Fisher #31985-070
ONE-Step™ Luciferase Assay System	BPS Bioscience #60690
Luminometer	

**Storage Conditions**

Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a -80°C freezer for long term storage. Contact technical support at support@bpsbioscience.com if the cells are not frozen in dry ice upon arrival.

**Media Formulations**

For best results, the use of validated and optimized media from BPS Bioscience is *highly recommended*. Other preparations or formulations of media may result in suboptimal performance.



Note: Thaw Media do *not* contain selective antibiotics. However, Growth Media *do* contain selective antibiotics, which are used to maintain selective pressure on the cell population expressing the gene of interest. Cells should be grown at 37°C with 5% CO<sub>2</sub>. BPS Bioscience's cell lines are stable for at least 10 passages when grown under proper conditions.

*Media Required for Cell Culture*

*Thaw Medium 1 (BPS Bioscience #60187):*

MEM medium supplemented with 10% FBS, 1% non-essential amino acids, 1 mM Na pyruvate, 1% Penicillin/Streptomycin.

*Growth Medium 1U (BPS Bioscience #78548):*

MEM medium supplemented with 10% FBS, 1% non-essential amino acids, 1 mM Na pyruvate, 1% Penicillin/Streptomycin, 0.5 µg/ml of Puromycin and 100 µg/ml of Hygromycin B.

*Media Required for Functional Cellular Assay*

*Assay Medium:*

Opti-MEM reduced serum medium.



GLP-1R agonists can be unstable in the presence of serum. Replace Thaw Medium 1 with Assay Medium if testing the agonist mentioned above. If the test agonist is stable in the presence of 10% Fetal Bovine Serum (FBS), Thaw Medium 1 may be used instead of Assay Medium.

**Cell Culture Protocol**

**Note: HEK293 cells are derived from human material and thus the use of adequate safety precautions is recommended.**

*Cell Thawing*

1. Retrieve a cell vial from liquid nitrogen storage. Keep on dry ice until ready to thaw.
2. When ready to thaw, swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. Once cells are thawed (it may be slightly faster or slower than 60 seconds), quickly transfer the entire content of the vial to an empty 50 ml conical tube.

**Note: Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.**

3. Using a 10 ml serological pipette, slowly add 10 ml of pre-warmed Thaw Medium 1 to the conical tube containing the cells. Thaw Medium 1 should be added dropwise while gently rocking the conical tube to permit gentle mixing and avoid osmotic shock.
4. Immediately spin down the cells at 300 x *g* for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed Thaw Medium 1.
5. Transfer the resuspended cells to a T25 flask or T75 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
6. After 24 hours of culture, check for cell attachment and viability. Change medium to fresh Thaw Medium 1 and continue growing in a 5% CO<sub>2</sub> incubator at 37°C until the cells are ready to passage.

*Note: Recovery of the frozen cells can take longer than a week. Change medium to fresh Thaw Medium 1 after a week.*

7. Cells should be passaged before they are fully confluent. At first passage and subsequent passages, use Growth Medium 1U.

### *Cell Passage*

1. Aspirate the medium, wash the cells with phosphate buffered saline (PBS) without  $\text{Ca}^{2+}/\text{Mg}^{2+}$ , and detach the cells from the culture vessel with 0.05% Trypsin/EDTA.
2. Once the cells have detached, add Growth Medium 1U and transfer to a tube.
3. Spin down the cells at  $300 \times g$  for 5 minutes, remove the medium and resuspend the cells in Growth Medium 1U.
4. Seed into new culture vessels at the recommended sub-cultivation ratio of 1:5 to 1:8 weekly or twice per week.

### *Cell Freezing*

1. Aspirate the medium, wash the cells with PBS without  $\text{Ca}^{2+}/\text{Mg}^{2+}$ , and detach the cells from the culture vessel with 0.05% Trypsin/EDTA.
2. Once the cells have detached, add Growth Medium 1U and count the cells.
3. Spin down the cells at  $300 \times g$  for 5 minutes, remove the medium and resuspend the cells in 4°C Cell Freezing Medium (BPS Bioscience #79796) at  $\sim 2 \times 10^6$  cells/ml.
4. Dispense 1 ml of cell suspension into each cryogenic vial. Place the vials in an insulated container for slow cooling and store at -80°C overnight.
5. Transfer the vials to liquid nitrogen the next day for long term storage.



Note: It is recommended to expand the cells and freeze at least 10 vials at an early passage for future use.

**Functional Validation**

- The following assays are designed for 96-well format. To perform the assay in a different format, the cell number and reagent volume should be scaled appropriately.
- The assay conditions should be performed in triplicate.
- The assay should include “Background Control”, “Stimulated”, and “Unstimulated Control” conditions.

*Assay Medium:* Opti-MEM reduced serum medium.

**A. Dose response of Rat GLP-1R/CRE Luciferase Reporter-HEK293 Cell Line to GLP-1R agonists**

1. Seed GLP-1R/CRE Luciferase Reporter HEK293 cells into a white clear-bottom 96-well cell culture plate at a density of ~30,000 cells per well in 90 µl of Assay Medium. Leave a few wells empty for use as the cell-free control wells (“Background Control”).
2. Incubate cells at 37°C in a CO<sub>2</sub> incubator for 16 to 24 hours.
3. The next day, prepare a serial dilution of GLP-1R agonists in Assay Medium at 10x testing concentration (10 µl/well). For a serial dilution, it is recommended to change a pipet tip for each dilution to avoid sample carry over.
4. Add 10 µl of the diluted GLP-1R agonist dilutions to the “Stimulated” wells.
5. Add 10 µl of Assay Medium to the “Unstimulated Control” wells.
6. Add 100 µl of Assay Medium to the “Background Control” wells (for determining background luminescence).
7. Incubate the plate at 37°C in a CO<sub>2</sub> incubator for ~ 5 hours.
8. Add 100 µl of the ONE-Step™ Luciferase reagent per well.
9. Rock gently at Room Temperature (RT) for ~15 minutes.
10. Measure luminescence using a luminometer.
11. Data Analysis: Subtract the average background luminescence (cell-free wells) from the luminescence reading of all wells. The fold induction of CRE luciferase reporter expression is the average background-subtracted luminescence of the stimulated wells divided by the average background-subtracted luminescence of unstimulated control wells.

*Note: Background subtracted luminescence of unstimulated cells is typically low in this cell line and can vary from assay to assay. This variability can lead to a broad range of fold induction values. Fold calculation can be calculated as follows:*

$$\text{Fold induction} = \frac{(\text{luminescence of stimulated cells} - \text{average background})}{(\text{average luminescence of unstimulated cells} - \text{average background})}$$

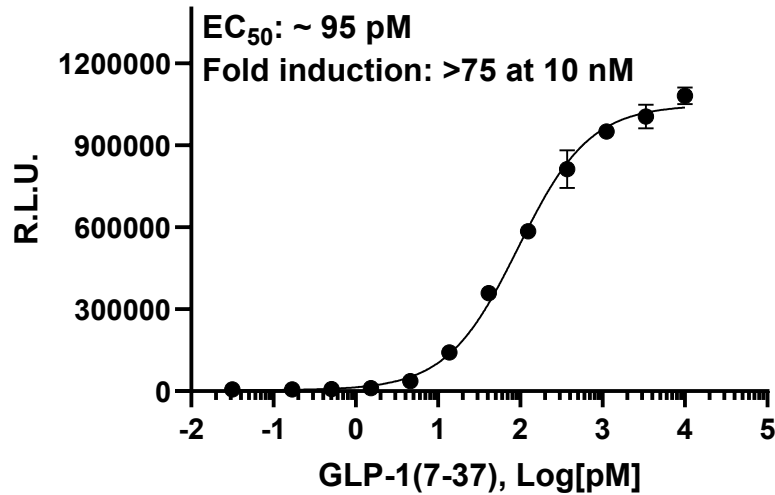


Figure 2. Dose response curve of Rat GLP-1R/CRE Luciferase Reporter HEK293 cells to GLP-1 (7-37). Cells were treated with increasing concentrations of GLP-1 (7-37). GLP-1 (7-37), which stimulated rat GLP-1R, inducing luciferase activity. Luciferase activity was measured with ONE-Step™ Luciferase Assay System. Results are expressed as relative luminescence versus the unstimulated control and the fold induction is calculated as described.

Data shown is representative.

### Sequence

Rattus Norvegicus GLP-1R sequence (NM\_012728.2)

MAVTPSLLRLALLLLGAVGRAGPRPQGATVSLSETVQKWREYRHQCQRFLTEAPLLATGLFCNRTFDDYACWPDGPPGSFVNVS  
CPWYLPWASSVLQGHVYRFCTAEGIWLHKDNSSLPWRDLSECEESKQGERNSPEEQLLSLYIIYTVGYALSFSALVIASAILVSFRHL  
HCTRNYIHLNLFASFILRALS VFIKDAALKWMYSTAAQQHQWDGLLSYQDSLGCRLVFLLMQYCVAANYWLLVEGVLYTLLAF  
SVFSEQRIFKLYLSIGWGVPLLFVIPWGVIVKLYEDEGCWTRNSNMNYWLIIRLPILFAIGVNFILFIRVICIVIAKLMCKTDIKC  
RLAKSTLTLIPLLGTHEVIFAFVMDEHARGTLRFVKLFTLSFTSQGFMVAVLYCFVNNEVQMEFRKSWERWRLERLNIQRDSS  
MKPLKCPTSSVSSGATVGSSVYAATCQNSCS

### References

Zhao X., et al., 2021 *Front. Endocrinol.* 12:721135.

### License Disclosure

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

### Troubleshooting Guide

Visit [bpsbioscience.com/cell-line-faq](https://bpsbioscience.com/cell-line-faq) for detailed troubleshooting instructions. For lot-specific information and all other questions, please visit <https://bpsbioscience.com/contact>.

**Related Products**

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
GLP-1R/CRE Luciferase Reporter HEK293 Cell Line	78176	2 vials
Rat GIPR/CRE Luciferase Reporter HEK293 Cell Line	83881	2 vials
GIPR/CRE Luciferase Reporter HEK293 Cell Line	78589	2 vials
Adenosine A2A Receptor Functional Cell Line	79381	2 vials
CGRPR/CRE Luciferase Reporter HEK293 Cell Line	78325	2 vials
CRE/CREB Luciferase Reporter HEK293 Cell Line	60515	2 vials

*Version 010926*