

DATA SHEET

CELL LINE DESIGNATION
ORIGIN (PARENTAL CELL)
GENE INTRODUCED
RECEPTOR INTRODUCED:

Glucagon-like peptide 1 receptor cell line (CB80200-248)
HEK 293-CNG cell (CB80200-200)
Genbank Locus ID 2740
Human glucagons-like peptide 1 receptor (NCBI protein database NP 002053.2)

USAGE

- cAMP assay for Gs-coupled human glucagons-like peptide 1 receptor.
- HEK293-CNG cells (CB80200-200) without transfected glucagons-like peptide 1 receptor are used as a negative control.

QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for glucagons-like peptide 1 receptor specific response.
3. Surviving rate: More than 2.5 million/vial on the second day after thawing.
4. The receptor specific activity is stable for 10 weeks continuous passage.

CELL CULTURE CONDITION

1. Growth medium: 90% DMEM, 10% FBS, 250 $\mu\text{g/ml}$ G418 and 1 $\mu\text{g/ml}$ puromycin
2. Freezing medium: 10% DMSO, 90% FBS

DATA EXAMPLE

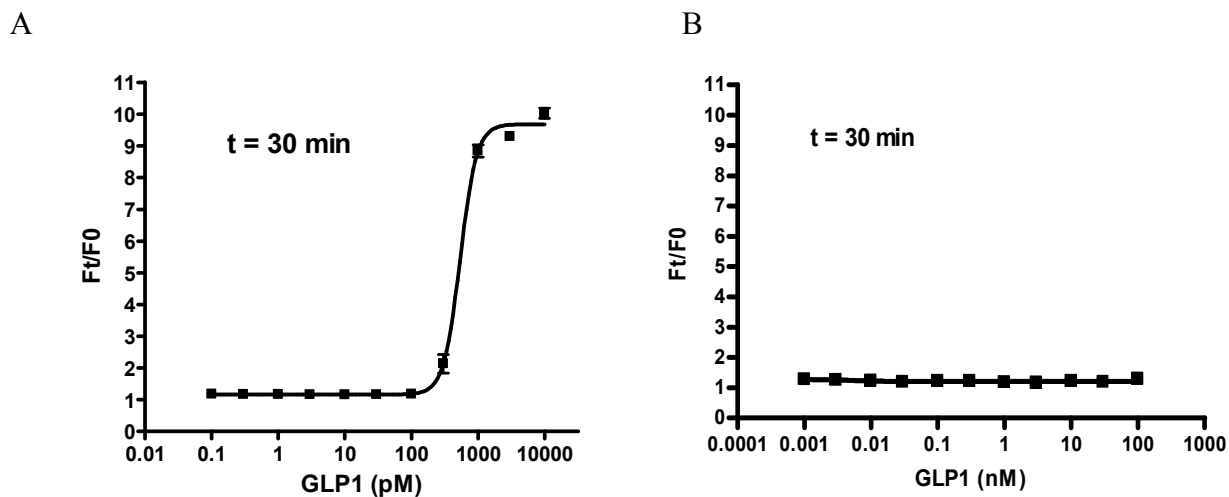


Figure 1. Response of ACTOne GLP1R cell line & parental cell line to GLP-1.

ACTOne GLP1R cells and parental cells (CB80200-200) were plated overnight in 20 μl culture medium on a BD Biocoat 384 well plate. The next day, cells were dye-loaded with 20 $\mu\text{l/well}$ of 1X Dye-loading solution (ACTOne Membrane Potential Assay Kit). After 2 hours of incubation at room temperature, two readings were obtained prior to and 30 min after the addition of GLP-1. Ratios of the two readings (F/F0) are plotted in the figure.

- A. Dose response curve of GLP-1 in ACTOne GLP1R cell line. EC50 = 536 pM in the presence of 25 μM of PDE inhibitor Ro20-1724.**
- B. Parental cells do not respond to GLP-1.**