

## DATA SHEET

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

Glucagon-like peptide 2 receptor cell line (CB-80200-241)  
HEK 293-CNG cell (CB-80200-200)  
Genbank Locus ID 9340  
Human glucagons-like peptide 2 receptor (NCBI protein database NP\_004237.1)

### USAGE

- cAMP assay for Gs-coupled human Glucagon-like peptide 2 receptor (GLP2R).
- HEK293-CNG cells (CB-80200-200) without transfected Glucagon-like peptide 2 receptor are used as a negative control.

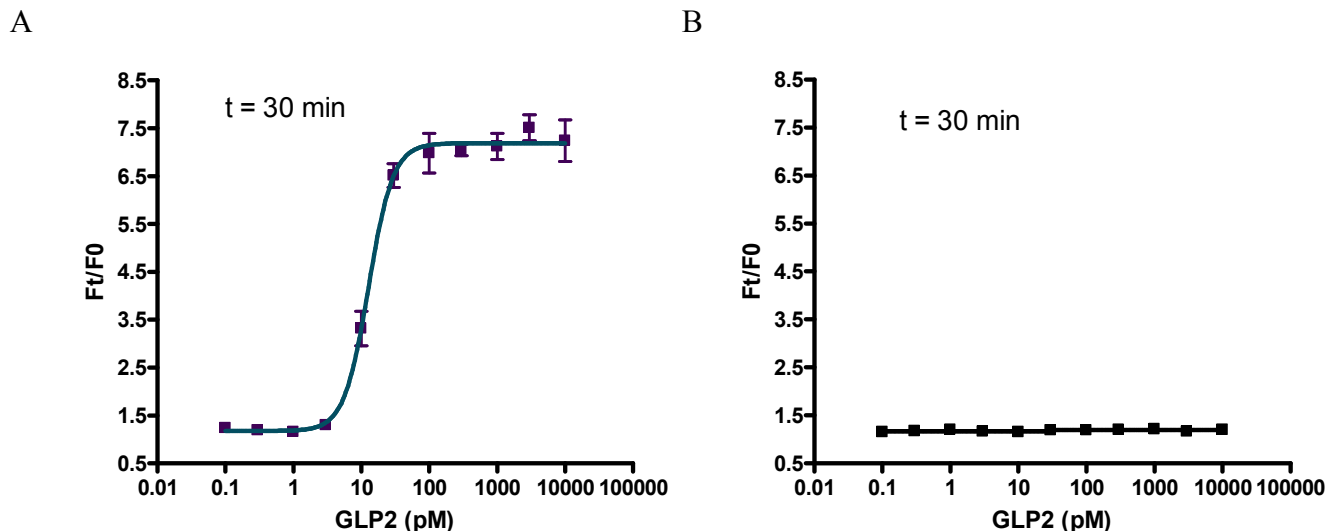
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Glucagon-like peptide 2 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 µg/ml G418 and 1 µg/ml puromycin
2. Freezing medium: 10% DMSO, 90% FBS

### DATA EXAMPLE



**Figure 1. Response of ACTOne GLP2R cell line & parental cell line to GLP-2.**

ACTOne GLP2R cells and parental cells (CB-80200-200) were plated overnight in 20 µl culture medium on a BD Biotac 384 well plate. The next day, cells were dye-loaded with 20 µl/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-2. Ratios of the two readings (Ft/F0) are plotted in the figure.

- A. Dose response curve of GLP-2 in ACTOne GLP2R cell line. EC50 = 12.8 pM in the presence of 25 µM of PDE inhibitor Ro20-1724.**
- B. Parental cells do not respond to GLP-2.**