

DATA SHEET

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

Prostaglandin E Receptor 4 cell line (CB-80200-219)
HEK 293-CNG cell (CB-80200-200)
Genbank Locus ID 5734
Human Prostaglandin E Receptor 4 (NCBI protein database NP_000949)

USAGE

- cAMP assay for Gs-coupled human Prostaglandin E Receptor 4 (PTGER4).
- HEK293-CNG cells (CB-80200-200) without transfected Prostaglandin E Receptor 4 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 Prostaglandin E Receptor 4 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

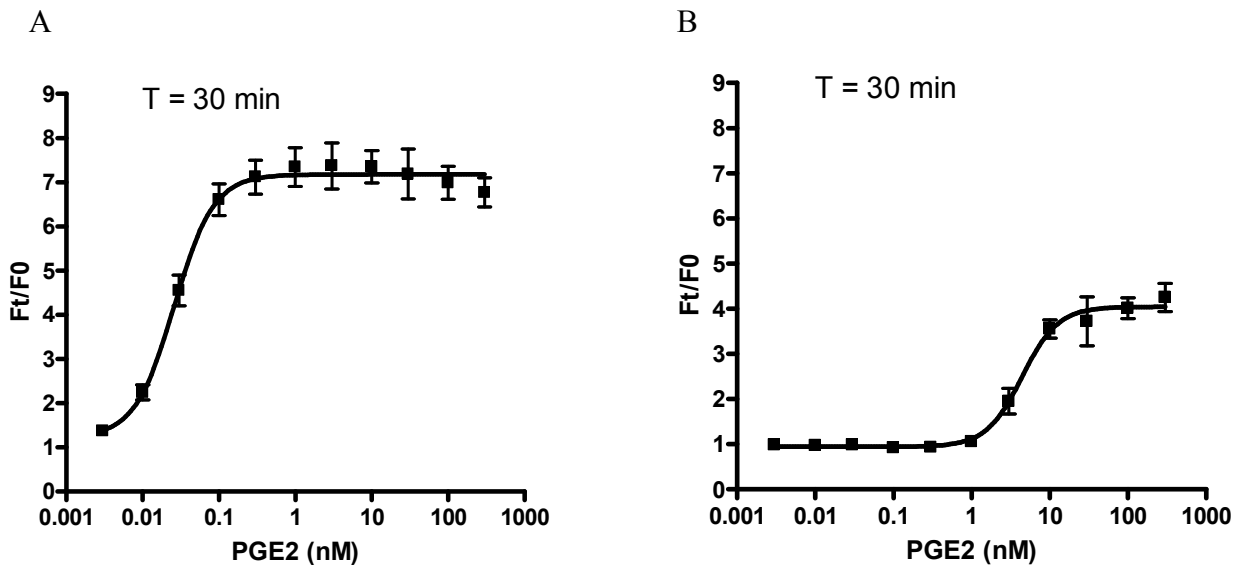


Figure 1. Response of ACTOne PTGER4 cell line & parental cell line to PGE2.

ACTOne PTGER4 cells and parental cells (CB-80200-200) were plated overnight in 20 µl culture medium on a BD Biocoat 384 well plate. The next day, cells were dye-loaded with 20 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 PGE2. Ratios of the two readings (F1/F0) are plotted in the figure.

- A. Dose response curve of PGE2 in ACTOne PTGER4 cell line. EC50 = 26.2 pM in the presence of PDE inhibitor Ro20-1724, and EC50 = 121 pM in the absence of Ro20-1724 (Data not shown).**
- B. Parental cells do not respond to PGE2 when the concentrations are less than 1 nM (in presence of PDE inhibitor). In the absence of PDE inhibitor, parental cells do not respond to PGE2 when the concentrations are less than 10 nM (Data not shown)**