

DATA SHEET

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

Prostaglandin I2 Receptor cell line (CB-80200-247)
HEK 293-CNG cell (CB-80200-200)
Genbank Locus ID 5739
Human Prostaglandin I2 Receptor (NCBI protein database
NP_000951.1)

USAGE

- cAMP assay for Gs-coupled human Prostaglandin I2 Receptor (PTGIR).
- HEK293-CNG cells (CB-80200-200) without transfected Prostaglandin I2 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 Prostaglandin I2 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

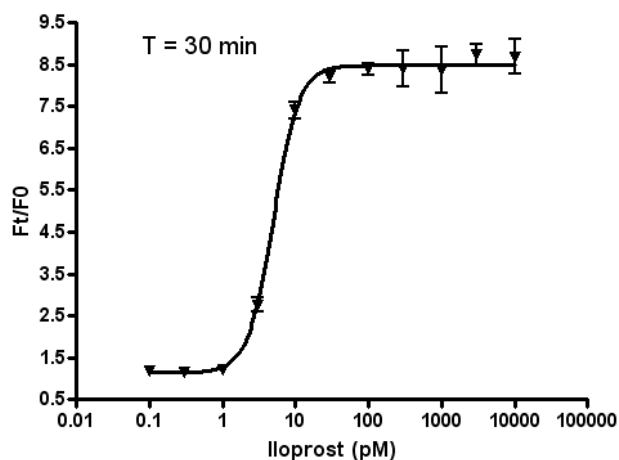


Figure 1. Response of ACTOne PTGIR cell line to Iloprost.

ACTOne PTGIR cells 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 Iloprost. Ratios of the two readings (Ft/F0) are plotted in the figure.

EC50 = 5.0 pM in the presence of PDE inhibitor Ro20-1724.

Parental cells don't respond to Iloprost (Data not shown)