

## DATA SHEET

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

Prostaglandin D Receptor 2 cell line (CB-80200-255)  
HEK 293-CNG cell (CB-80200-200)  
Genbank Locus ID 5729  
Human Prostaglandin D Receptor 2 (NCBI protein database NP\_000944.1)

### USAGE

- cAMP assay for Gs-coupled human Prostaglandin D Receptor 2 (PTGDR2).
- HEK293-CNG cells (CB-80200-200) without transfected Prostaglandin D Receptor 2 are used as a negative control.

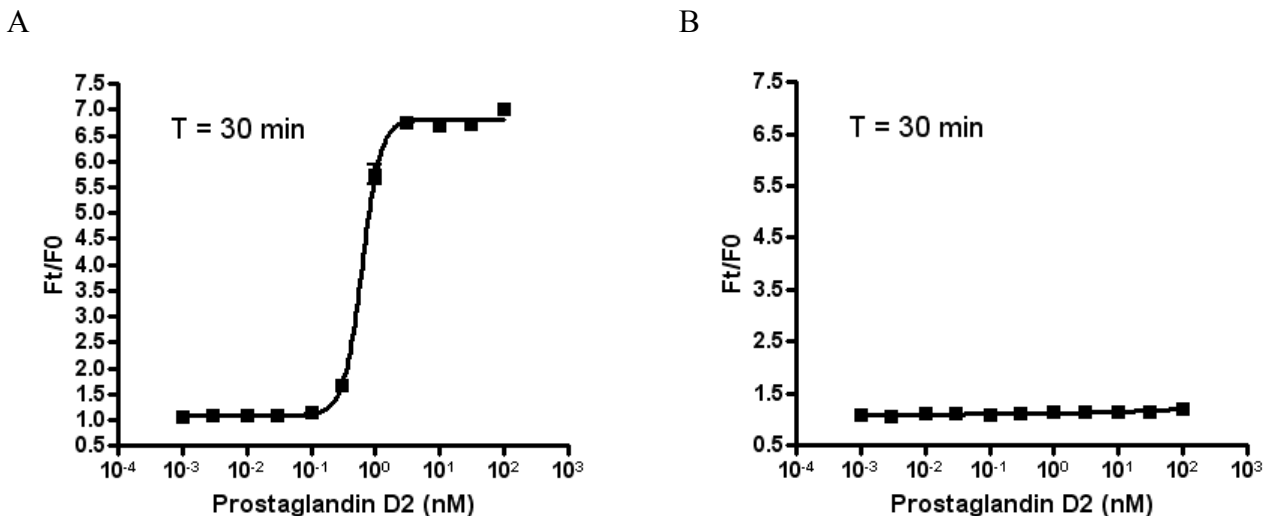
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Prostaglandin D Receptor 2 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 Prostaglandin D Receptor 2 cell line & parental cell line to prostaglandin D2.**

ACTOne PTGDR2 cells and parental cells (CB-80200-200) were plated overnight in 20  $\mu\text{l}$  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 prostaglandin D2. Ratios of the two readings (F/F<sub>0</sub>) are plotted in the figure.

- A. Dose response curve of prostaglandin D2 in ACTOne PTGDR2 cell line. EC<sub>50</sub> = 0.62 nM in the presence of PDE inhibitor Ro20-1724.**
- B. Parental cells do not respond to prostaglandin D2.**