

DATA SHEET

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

Phosphodiesterase 1B (cAMP) cell line (CB-81200-103)
HEK 293-CNG-Gs cell (CB-81200-100)
Genbank Locus ID 5153
Human phosphodiesterase 1B (NCBI protein database NP_000915.1)

USAGE

- cAMP assay for recombinant human phosphodiesterase 1B (PDE1B).
- HEK293-CNG-Gs cells (CB-81200-100) 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 PDE1B specific response.
3. Surviving rate: More than 2.5 million/vial on the second day after thawing.

CELL CULTURE CONDITION

1. DMEM-10%FBS supplemented with 250 µg/ml G418, 1 µg/ml Puromycin and 5 µg/ml blasticidin.
2. Freezing medium: 10% DMSO, 90% complete cell culture medium.

DATA EXAMPLE

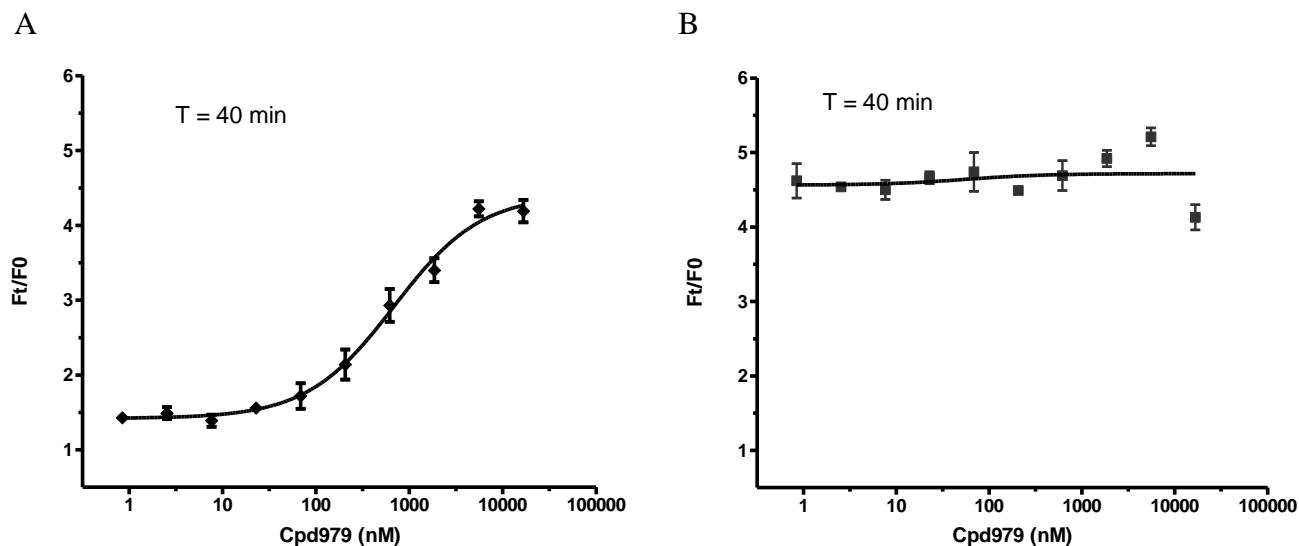


Figure 1. Response of ACTOne PDE1B cell line & parental cell line to Cpd979.

ACTOne PDE1B cells and parental cells (CB-81200-100) were plated overnight in 20 µl culture medium on a 384 well Biocoat plate. The next day, cells were dye-loaded with 20 µl/well of ACTOne membrane potential dye. After 2 hour of incubation at room temperature, baseline was recorded using a FlexStation (Molecular Devices) (F0). 10 µl of PDE inhibitors at various concentrations (with 0.5 µM Forskolin) were added to the cell plate, and the data was recorded 40 minutes (Ft) after drug addition. Dose response curves were generated by Prism.

- A. Dose response curve of Cpd979 in ACTOne PDE1B cell line. IC50 = 684 nM in the presence of 0.1 µM Forskolin**
- B. Parental cells do not respond to Cpd979 in the presence of 0.1 µM Forskolin**