

DATA SHEET

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

Phosphodiesterase 2A cell line (CB-81200-101)
HEK 293-CNG-Gs cell (CB-81200-100)
Genbank Locus ID 5138
Human phosphodiesterase 2A (NCBI protein database NP_001137311.1)

USAGE

- cAMP assay for recombinant human phosphodiesterase 2A (PDE2A).
- 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 PDE2A 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. 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.
- 3.

DATA EXAMPLE

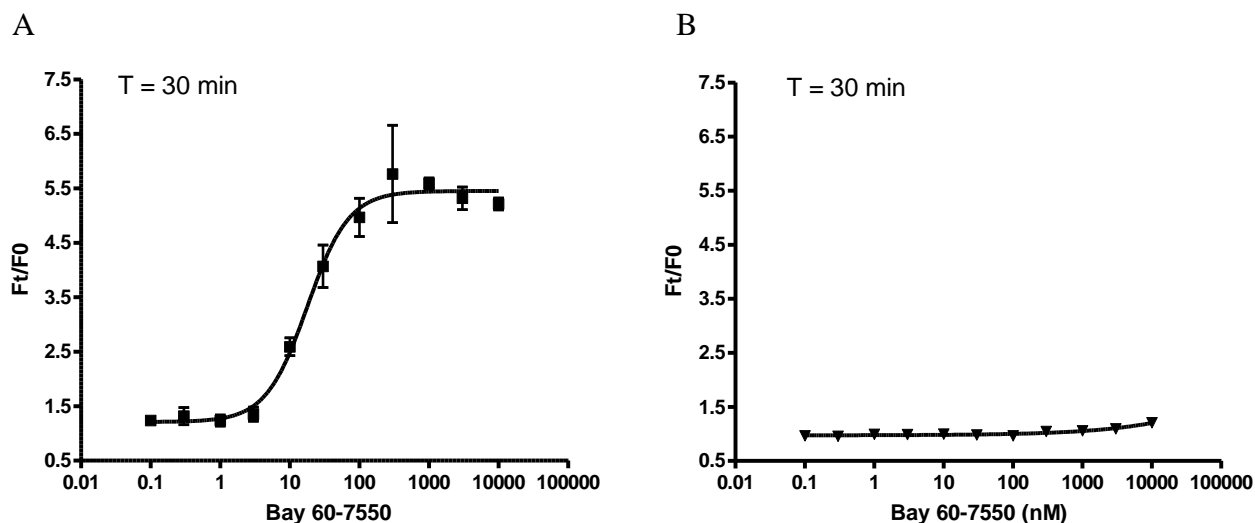


Figure 1. Response of ACTOne PDE2A cell line & parental cell line to Bay 60-7550.

ACTOne PDE2A 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 were added to the cell plate, and the data was recorded 30 minutes (Ft) after drug addition. Dose response curves were generated by Prism.

- A. Dose response curve of Bay 60-7550 in ACTOne PDE2A cell line. IC50 = 17.8 nM in the presence of 10 µM of Ro20-1724**
- B. Parental cells do not respond to Bay 60-7550 in the presence of 10 µM of Ro20-1724**