

## DATA SHEET

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

Phosphodiesterase 10A cell line (CB-81200-102)  
HEK 293-CNG-Gs cell (CB-81200-100)  
Genbank Locus ID 10846  
Human phosphodiesterase 10A (NCBI protein database  
AAD32595.1)

### USAGE

- cAMP assay for recombinant human phosphodiesterase 10A (PDE10A).
- HEK293-CNG-Gs cells (CB-81200-100) are used as a negative control.

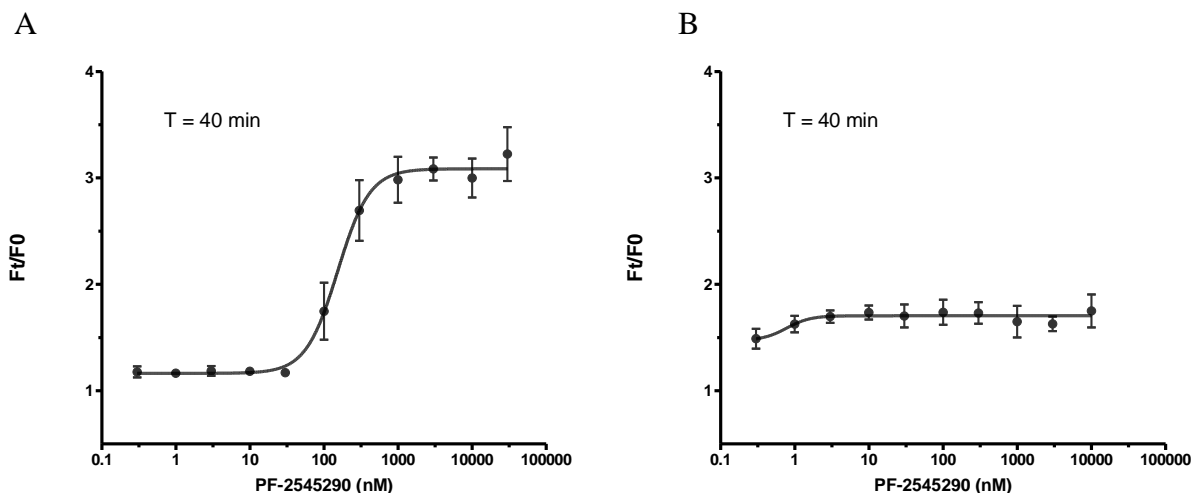
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for PDE10A 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.

### DATA EXAMPLE



**Figure 1. Response of ACTOne PDE10A cell line & parental cell line to PF-2545290.**

ACTOne PDE10A 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 40 minutes (Ft) after drug addition. Dose response curves were generated by Prism.

- A. Dose response curve of PF-2545290 in ACTOne PDE10A cell line. IC50 = 153 nM in the presence of 10 µM of Ro20-1724**
- B. Parental cells do not respond to PF-2545290 in the presence of 10 µM of Ro20-1724**