

APPENDIX

CELL LINE DESIGNATION

G protein-coupled receptor 119 cell line
(CB-80200-274)

ORIGIN (PARENTAL CELL)

HEK 293-CNG cell (CB80200-200)

GENE INTRODUCED

Genbank Locus ID 139760

RECEPTOR INTRODUCED:

Human G protein-coupled receptor 119
(NCBI protein database XP_066873.1)

USAGE

- cAMP assay for Gs-coupled human G protein-coupled receptor 119 .
- HEK293-CNG cells (CB80200-200) without transfected G protein-coupled receptor 119 is used as a negative control.

QUALITY CONTROL

1. This cell line has been tested positive for G protein-coupled receptor 119 specific response.
2. Surviving rate: More than 2.5 million/vial on the second day after thawing.

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% growth medium.

DATA EXAMPLE

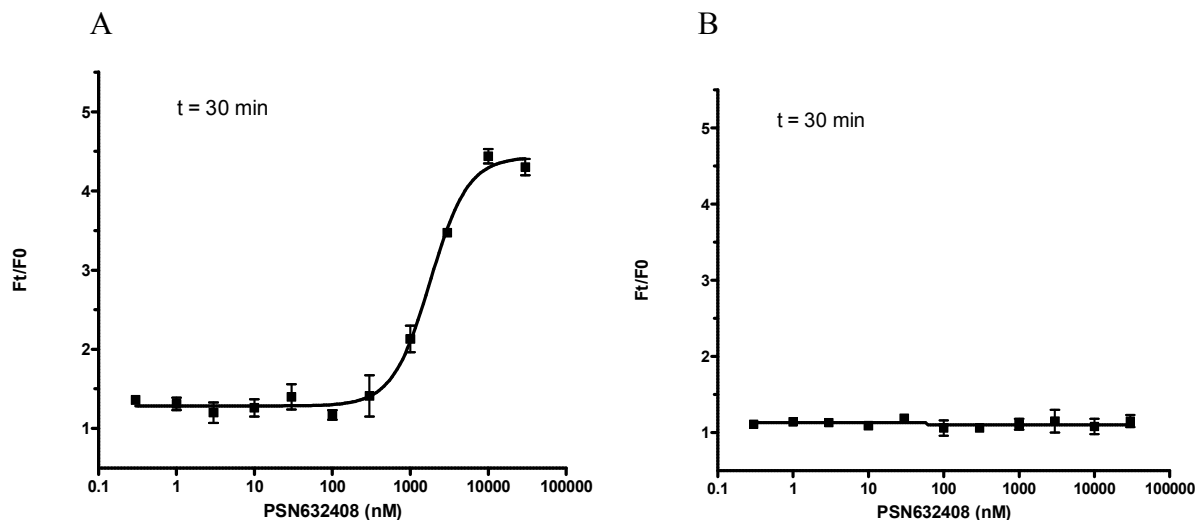


Figure 1. Response of ACTOne GPR119 cell line & parental cell line to PSN632408.

ACTOne GPR119 cells and parental cells (CB80200-200) 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 PSN632408. Ratios of the two readings (F/F0) are plotted in the figure.

- A. Dose response curve of PSN632408 in ACTOne GPR119 cell line. EC50 = 1.8 μM in the presence of 25 μM of PDE inhibitor Ro20-1724.**
- B. Parental cells do not respond to PSN632408.**