

DATA SHEET

CELL LINE DESIGNATION

Endothelial Differentiation, Sphingolipid G-Protein-coupled Receptor, 1 cell line (CB-80300-254)

ORIGIN (PARENTAL CELL)

HEK 293-CNG cell (CB-80200-200)

GENE INTRODUCED

Genbank Locus ID 1901

RECEPTOR INTRODUCED:

Human endothelial differentiation, sphingolipid G-protein-coupled receptor, 1 (NCBI protein database NP_001391.2)

USAGE

- cAMP assay for Gi-coupled human Endothelial Differentiation, Sphingolipid G-Protein-coupled Receptor, 1 (EDG1)
- HEK293-CNG cells (CB-80200-200) without transfected Endothelial Differentiation, Sphingolipid G-Protein-coupled Receptor, 1 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 Endothelial Differentiation, Sphingolipid G-Protein-coupled Receptor, 1 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 µg/ml G418 and 1 µg/ml puromycin
2. Freezing medium: 10% DMSO, 90% FBS

DATA EXAMPLE

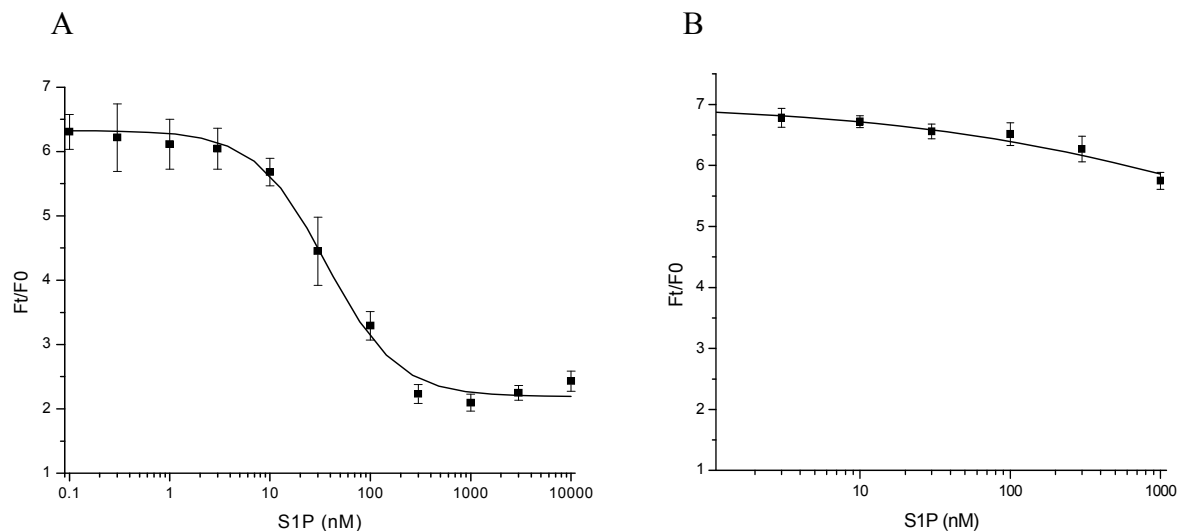


Figure 1. Response of ACTOne EDG1 cell line & parental cell line to S1P.

ACTOne EDG1 cells and parental cells (CB-80200-200) were plated overnight in 20 µl culture medium on a BD Biotec 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 40 min after the addition of S1P. Ratios of the two readings (F/F0) are plotted in the figure.

- A. Dose response curve of S1P in ACTOne EDG1 cell line. EC50 = 36 nM in the presence of PDE inhibitor Ro20-1724 and β-adrenoceptor agonist isoproterenol.**
- B. Parental cells do not respond to S1P.**