

DATA SHEET

CELL LINE DESIGNATION

Chemokine (C-X-C motif) Receptor 4 cell line
(CB-80300-265)

ORIGIN (PARENTAL CELL)

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

GENE INTRODUCED

Genbank Locus ID 7852

RECEPTOR INTRODUCED:

Human chemokine (C-X-C motif) receptor 4 (NCBI protein
database AF052572)

USAGE

- cAMP assay for Gi-coupled human Chemokine (C-X-C motif) Receptor 4 (CXCR4).
- HEK293-CNG cells (CB-80200-200) without transfected Chemokine (C-X-C motif) Receptor 4 are used as a negative control.

QUALITY CONTROL

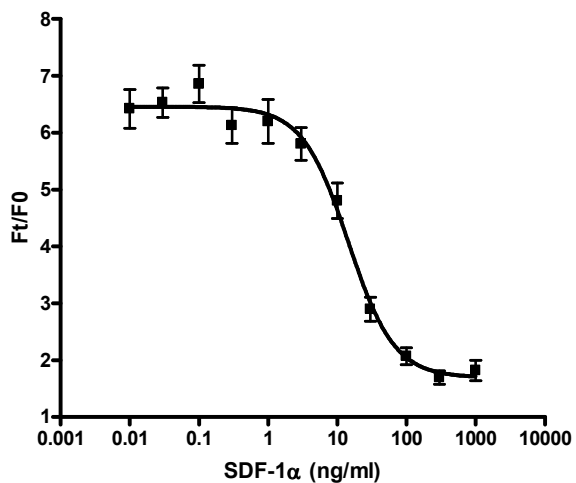
1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Chemokine (C-X-C motif) Receptor 4 specific response.
3. Surviving rate: More than 2.5 million/vial on the second day after thawing.
4. The receptor specific activity is stable for 10 weeks 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

A



B

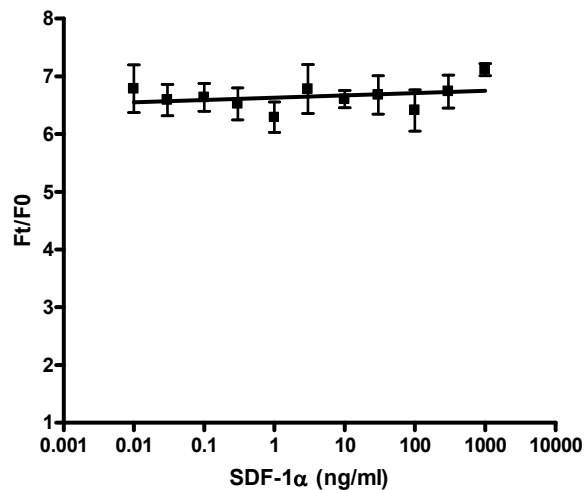


Figure 1. Response of ACTOne CXCR4 cell line & parental cell line to SDF-1α.

ACTOne CXCR4 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 25 min after the addition of SDF-1α. Ratios of the two readings (F/F0) are plotted in the figure.

- A. Dose response curve of SDF-1α in ACTOne CXCR4 cell line. EC50 = 14 ng/ml in the presence of PDE inhibitor Ro20-1724 and β-adrenoceptor agonist isoproterenol.**
- B. Parental cells do not respond to SDF-1α.**