

DATA SHEET

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

Arginine Vasopressin Receptor 2 cell line (CB-80200-246)
HEK 293-CNG cell (CB-80200-200)
Genbank Locus ID 554
Human Arginine Vasopressin Receptor 2 (NCBI protein database NP_000045.1)

USAGE

- cAMP assay for Gs-coupled human Arginine Vasopressin Receptor 2 (AVPR2).
- HEK293-CNG cells (CB-80200-200) without transfected Arginine Vasopressin Receptor 2 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 Arginine Vasopressin Receptor 2 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

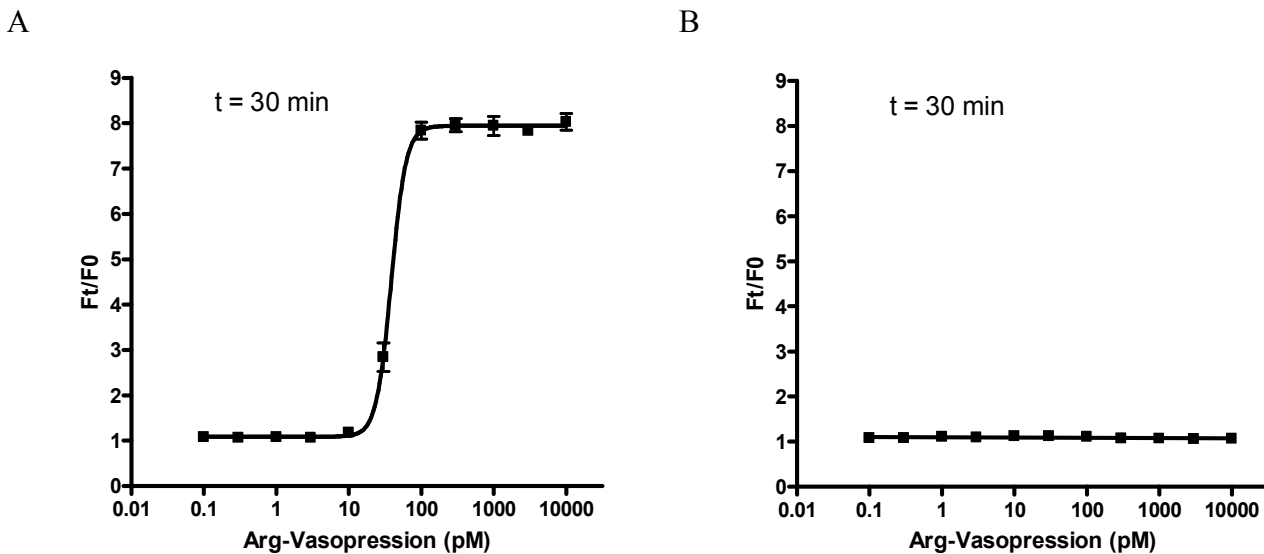


Figure 1. Response of ACTOne AVPR2 cell line & parental cell line to Arginine Vasopressin.

ACTOne AVPR2 cells and parental cells (CB-80200-200) were plated overnight in 20 µl culture medium on a BD Biotac 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 Arginine Vasopressin. Ratios of the two readings (F_t/F_0) are plotted in the figure.

- A. Dose response curve of Arginine Vasopressin in ACTOne AVPR2 cell line. $EC_{50} = 38.6$ pM in the presence of PDE inhibitor Ro20-1724, and $EC_{50} = 110$ pM in the absence of Ro20-1724 (data not shown).**
- B. Parental cells do not respond to Arginine Vasopressin.**