

## DATA SHEET

### CELL LINE DESIGNATION

Adenylate Cyclase Activating Polypeptide 1 Receptor cell line (CB-80200-234)

### ORIGIN (PARENTAL CELL)

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

### GENE INTRODUCED

Genbank Locus ID 117

### RECEPTOR INTRODUCED:

Human Adenylate Cyclase Activating Polypeptide 1 Receptor (NCBI protein database P41586)

### USAGE

- cAMP assay for Gs-coupled human Adenylate Cyclase Activating Polypeptide 1 Receptor (ADCYAP1R1).
- HEK293-CNG cells (CB-80200-200) without transfected Adenylate Cyclase Activating Polypeptide 1 Receptor are used as a negative control.

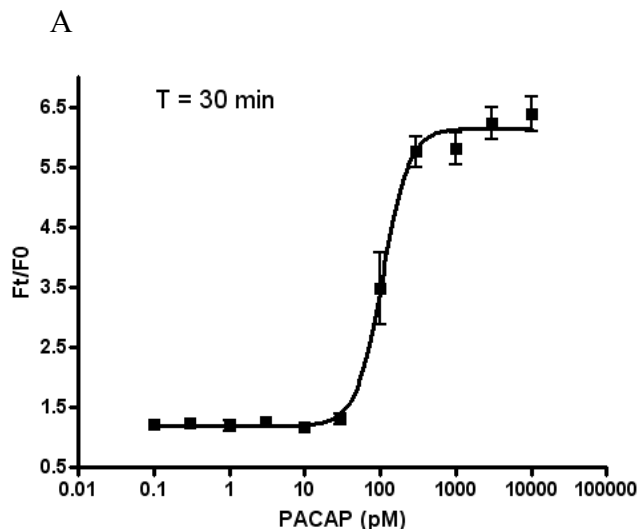
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Adenylate Cyclase Activating Polypeptide 1 Receptor 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 ADCYAP1R1 cell line to PACAP.

ACTOne ADCYAP1R1 cells 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 PACAP. Ratios of the two readings (Ft/F0) are plotted in the figure.

- A. Dose response curve of PACAP in ACTOne ADCYAP1R1L cell line. EC50 = 107.5 pM in the absence of PDE inhibitor Ro20-1724.
- B. Parental cells do not respond to PACAP (data not shown).