

DATA SHEET

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

Adenosine A2b receptor cell line (CB-80200-260)
HEK 293-CNG cell (CB-80200-200)
Genbank Locus ID 136
Human Adenosine A2b receptor (NCBI protein database
NP_000667.1)

USAGE

- cAMP assay for Gs-coupled human Adenosine A2b receptor (ADORA2B).
- HEK293-CNG cells (CB-80200-200) without transfected Adenosine A2b 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 Adenosine A2b 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

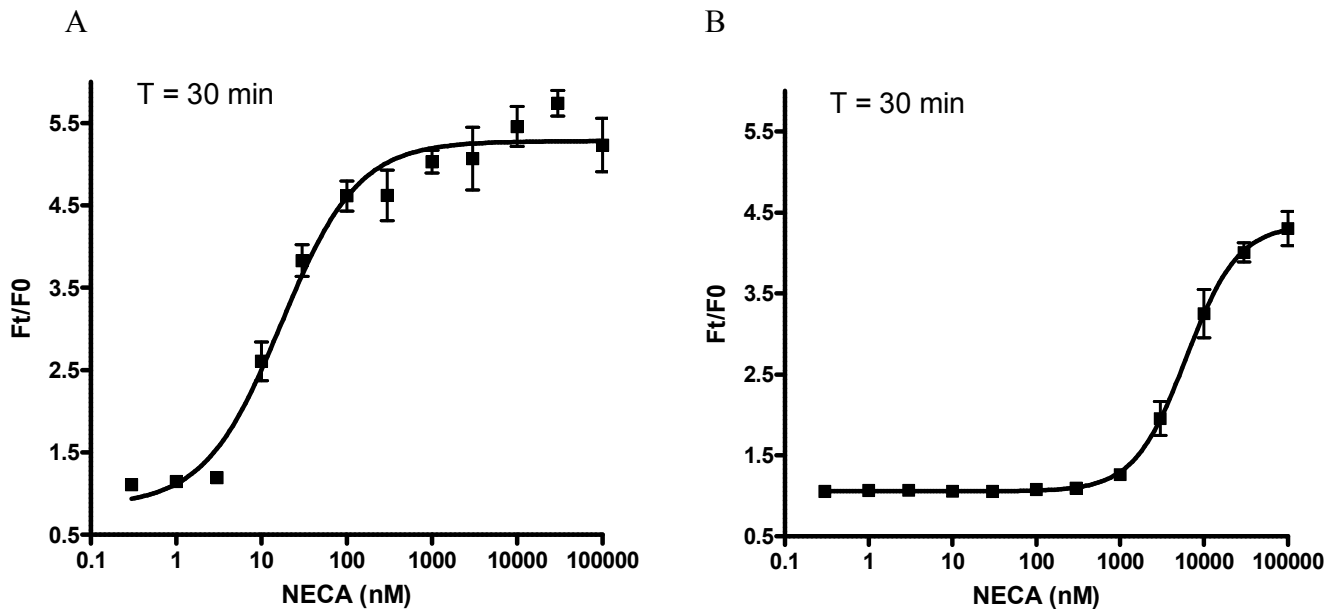


Figure 1. Response of ACTOne ADORA2B cell line & parental cell line to NECA.

ACTOne ADORA2B 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 30 min after the addition of NECA. Ratios of the two readings (Ft/F0) are plotted in the figure.

- A. Dose response curve of NECA in ACTOne ADORA2B cell line. EC50 = 17.08 nM in the absence of PDE inhibitor Ro20-1724.**
- B. Dose response curve of NECA in the parental cell line. EC50 = 6.1 µM in the absence of PDE inhibitor Ro20-1724.**