

## DATA SHEET

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

Adenosine A1 Receptor cell line (CB-80300-212)  
HEK 293-CNG cell (CB-80200-200)  
Genbank Locus ID 134  
Human Adenosine A1 Receptor (NCBI protein database  
NP\_000665)

### USAGE

- cAMP assay for Gi-coupled human Adenosine A1 Receptor (ADORA1).
- HEK293-CNG cells (CB-80200-200) without transfected Adenosine A1 Receptor are used as a negative control.

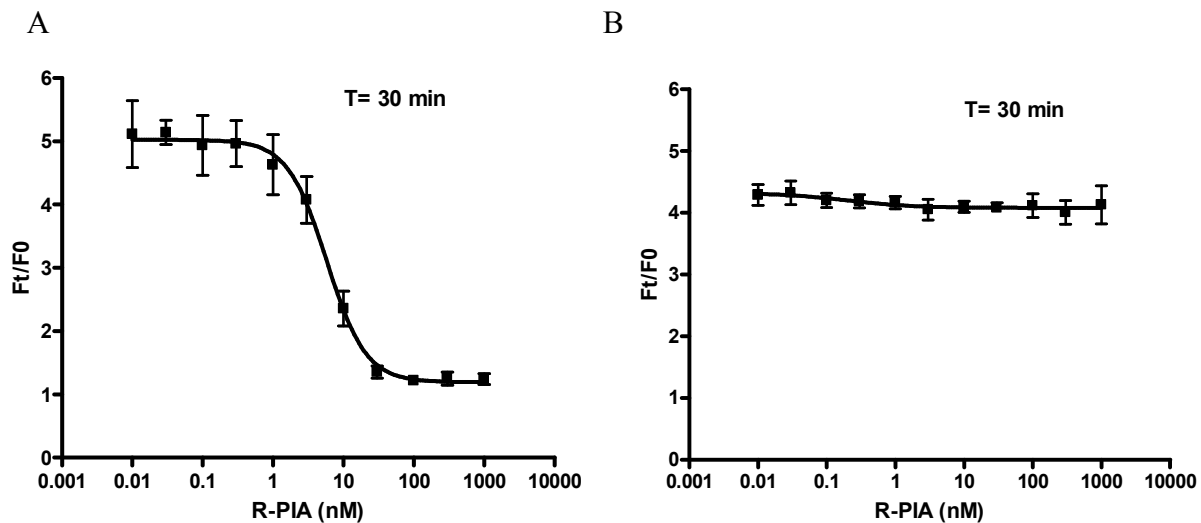
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Adenosine A1 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 ADORA1 cell line & parental cell line to R-PIA.

ACTOne ADORA1 cells and parental cells (CB-80200-200) 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 R-PIA. Ratios of the two readings (F/F<sub>0</sub>) are plotted in the figure.

- A. Dose response curve of R-PIA in ACTOne ADORA1 cell line. EC<sub>50</sub> = 5.8 nM in the presence of PDE inhibitor Ro20-1724 and β-adrenoceptor agonist isoproterenol.**
- B. Parental cells do not respond to R-PIA.**