

## DATA SHEET

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

Amylin 3 Receptor cell line (CB-80200-271)  
ACTOne Calcitonin Receptor cells (CB-80200-258)  
EMBL Accession AJ001016.1  
RAMP3 (NCBI protein database CAA04474)

### USAGE

- cAMP assay for Gs-coupled human Amylin 3 Receptor (AMY3).
- ACTOne Calcitonin Receptor cells (CB-80200-258) without transfected RAMP3 are used as a negative control.

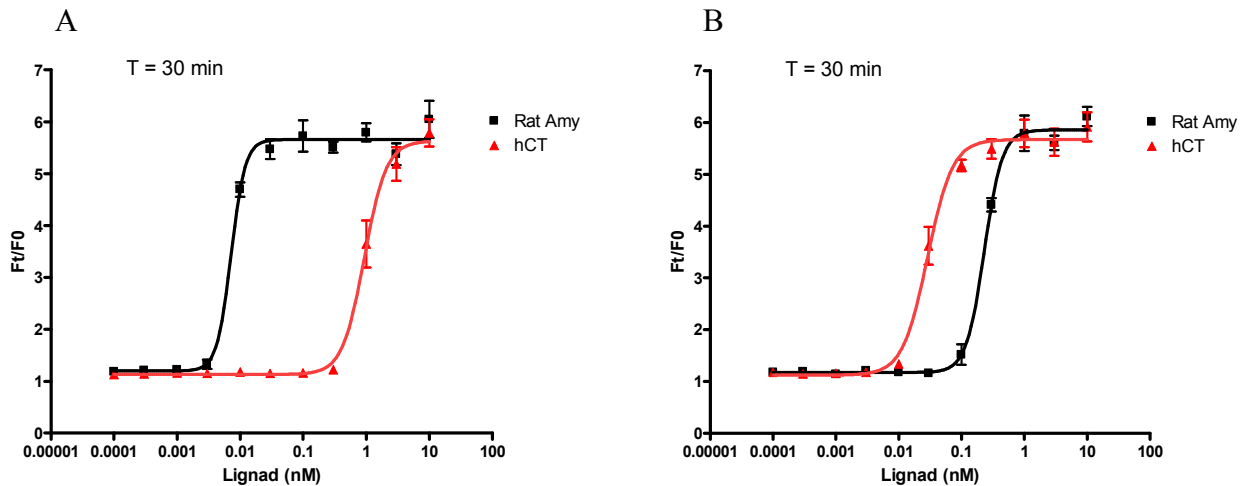
### QUALITY CONTROL

1. This cell line has been tested negative for *Mycoplasma sp.*
2. This cell line has been tested positive for Amylin 3 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, 1 µg/ml puromycin and 150 µg/ml hygromycin B
2. Freezing medium: 10% DMSO, 90% FBS

### DATA EXAMPLE



#### Figure 1. Response of ACTOne AMY3 cell line & parental cell line to amylin and calcitonin.

ACTOne AMY3 cells and parental cells (CB-80200-258) 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 rat amylin or hCT. Ratios of the two readings (F/F0) are plotted in the figure.

- A. Dose response curve of rat amylin and human calcitonin in ACTOne AMY3 cell line. In the presence of PDE inhibitor Ro20-1724, EC50 = 7.2 pM with rAmy and EC50 = 934 pM with hCT.**
- B. Dose response curve of rat amylin and human calcitonin in Parental cells. In the presence of PDE inhibitor Ro20-1724, EC50 = 230 pM with rAmy and EC50 = 28.7 pM with hCT**