

DATA SHEET

CELL LINE DESIGNATION

5-Hydroxytryptamine (Serotonin) Receptor 4 cell line
(CB-80200-231)

ORIGIN (PARENTAL CELL)

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

GENE INTRODUCED

Genbank Locus ID 3360

RECEPTOR INTRODUCED:

Human 5-hydroxytryptamine (serotonin) receptor 4 (NCBI
protein database NP_000861.1)

USAGE

- cAMP assay for Gs-coupled human 5-Hydroxytryptamine (Serotonin) Receptor 4 (HTR4).
- HEK293-CNG cells (CB-80200-200) without transfected 5-Hydroxytryptamine (Serotonin) Receptor 4 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 5-Hydroxytryptamine (Serotonin) Receptor 4 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. Growth medium for ACTOne Assay: 90% DMEM, 10% FBS (CHARCOAL STRIPPED), 250 µg/ml G418 and 1 µg/ml puromycin
3. Freezing medium: 10% DMSO, 90% FBS