

DATA SHEET

CELL LINE DESIGNATION

Human Thyrotropin-releasing Hormone Receptor cell line (CB-80400-285)

ORIGIN (PARENTAL CELL)

HEK293

GENE INTRODUCED

Genbank Locus ID 7201

RECEPTOR INTRODUCED:

Human Thyrotropin-releasing Hormone Receptor (NCBI protein database NP_003292.1)

USAGE

- Calcium influx assay for Gq-coupled human Thyrotropin-releasing Hormone Receptor (TRHR).
- HEK293 cells without transfected Thyrotropin-releasing Hormone 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 Thyrotropin-releasing Hormone Receptor specific response.
3. Surviving rate: More than 2.5 million/vial on the second day after thawing.

CELL CULTURE CONDITION

1. Growth medium: 90% DMEM, 10% FBS and 1 µg/ml puromycin
2. Freezing medium: 10% DMSO, 90% FBS

DATA EXAMPLE

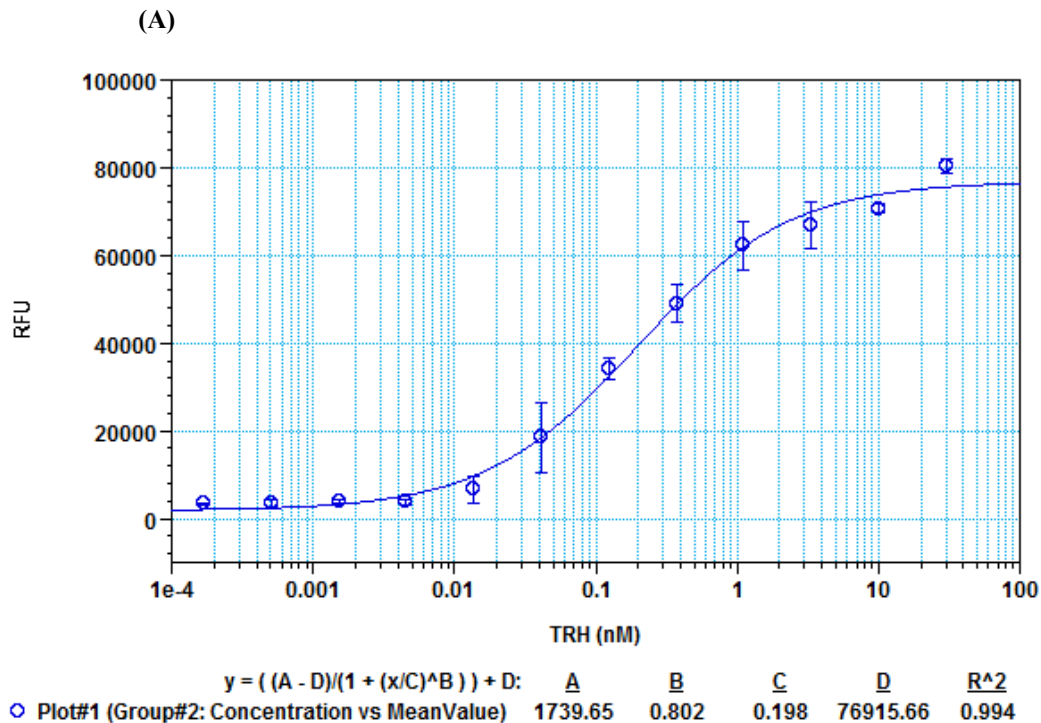


Figure 1. Response of HEK293-TRHR cell line to TRH.

TRHR cells and parental cells (HEK293) were plated overnight in 80 µl culture medium on a 96-well plate black/clear plate. The next day, the cells were loaded with equal volume of Codex non-wash calcium dye (CB-80500-301) and incubated at a 37°C incubator for 60 min. The plate was taken out and kept at room temperature for another 30 min. The assay was performed on a FlexStation with online addition of different concentrations of TRH

- A. Dose response curve of TRH in HEK293-TRH cell line. EC50 = 0.198 nM**
- B. Parental cells do not respond to TRH (data not shown).**