

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CT RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1231-1B

**Lot Number:** C1231-1B-031318

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Cellbanker 2 (Amsbio #11891)

**Host cell:** CHO-K1

**Transfection:** Full-length Human CALCR cDNA (GenBank Accession Number NM\_001742) with FLAG-tag sequence at the N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

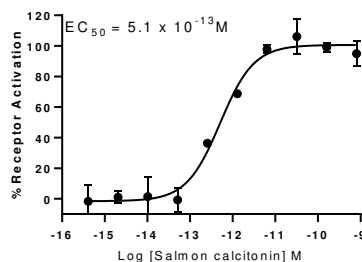
**Propagation Medium:** DMEM/F-12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** Stable for a minimum of two months in continuous culture

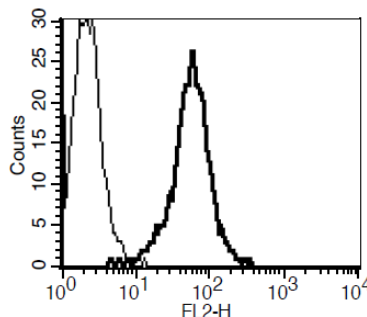
**Background:** Calcitonin receptor (CT or CALCR) is a receptor for calcitonin. In the presence of the receptor activity modifying proteins (RAMP1, 2 or 3), it acts as a receptor for amylin. CT is expressed in brain, bone, stomach, intestine, kidney and testes. Data from transgenic mice support a major role for calcitonin and its receptor in the formation and resorption aspects of bone metabolism under physiological conditions.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent increase of intracellular cAMP level upon treatment with ligand with transiently transfected cells, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

- Gorn *et al.* (1992) Cloning, characterization, and expression of a human calcitonin receptor from an ovarian carcinoma cell line. *J Clin Invest* 90:1726-1735.
- Taboulet *et al.* (1998) Calcitonin receptor polymorphism is associated with a decreased fracture risk in post-menopausal women. *Hum Mol Genet* 7:2129-2133.

**FOR RESEARCH USE ONLY.**

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.