

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CT RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1231-1B

Lot Number: C1231-1B-031318

Quantity: 1 vial (2×10^6) 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, 10% FBS, 10 μ g/mL puromycin

Stability: In progress

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

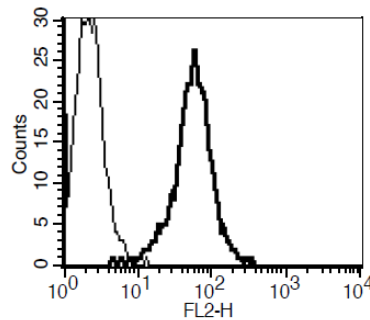


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand with transiently transfected cells, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG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