

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT CT RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MH1231

Lot Number: MH1231-021618

Quantity: 1 vial (4.4mg/mL, 1mg)

Packaging Buffer: 20mM Gly-Gly,
1mM MgCl₂, 25mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Full-length Human
CALCR cDNA (GenBank Accession
Number NM_001742) with FLAG-tag
sequence at the N-terminus

Recommended Storage: Store at
-80°C. Avoid repeated freeze-thaw.

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: Binding assays

Figure 1

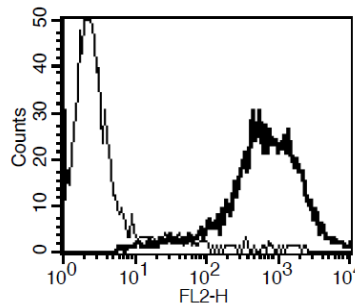


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.