

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT CASR RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1233

Lot Number: 08/13/10

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human CASR cDNA (GenBank accession number NM_000388) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: CASR is a calcium-sensing receptor and plays an important role in regulating PTH secretion. It is expressed in many different tissues, such as parathyroid cells, pituitary cells, kidney, fibroblasts, keratinocytes and human colon epithelial cells. CASR is a potential therapeutic target for the treatment of many diseases, including hyperparathyroidism and osteoporosis. Mutations in the CASR gene can result in gain or loss of receptor function. Familial Hypocalcemic Hypercalcemia (FHH) and Neonatal Severe Primary Hyperparathyroidism (NSHPT) have been associated with loss of CASR function, while Autosomal Dominant Hypocalcemia (ADH) and Bartter syndrome type V have been associated with gain of CASR function.

Application: Functional assays

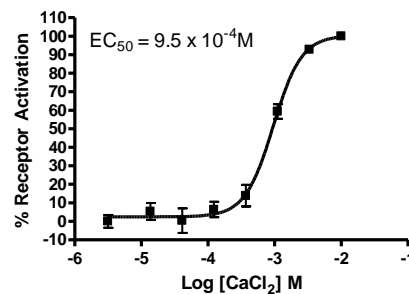


Figure legend: Dose-dependent accumulation of intracellular IP1 upon treatment with ligand, measured with IP-one Tb kit (Cisbio 62IPAPEC).

References:

D'Souza-Li (2006) The calcium-sensing receptor and related diseases. *Arq Bras Endocrinol Metabol* 50:628-639.

Romoli *et al.* (1999) Expression of calcium-sensing receptor and characterization of intracellular signaling in human pituitary adenomas. *J Clin Endocrinol Metab* 84:2848-2853.

FOR RESEARCH USE ONLY.

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