

## MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT AMY1 RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** DC1509-1a

**Lot Number:** DC1509-1a-040925

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2

**Host cell:** CHO-K1

**Transfection:** Full-length Human CALCR cDNA (GenBank Accession Number NM\_001742) with FLAG-tag sequence at the N-terminus and Full-length Human receptor activity modifying protein 1 (RAMP1) cDNA (GenBank Accession Number BC000548) with myc-tag at the C-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

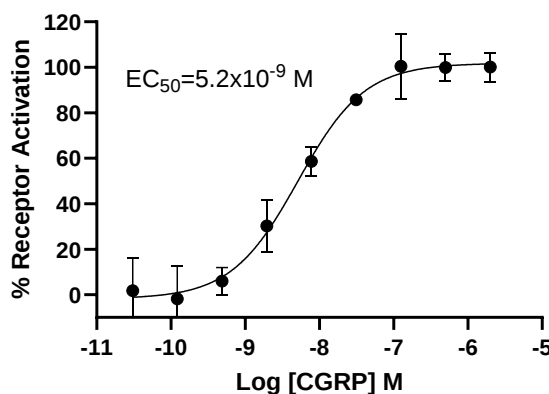
**Propagation Medium:** DMEM/F12, 10% FBS

### Data sheet

**Background:** Amylin receptors are multimeric complexes and have been closely associated with Calcitonin receptor (CT). CT, in the presence of the receptor activity modifying proteins (RAMP1, RAMP2 and RAMP3), forms pharmacologically distinct Amylin subtypes (AMY1, AMY2 and AMY3) which acts as a high affinity receptor for amylin, a hormone secreted by B cell of pancreas that has a major role in glucose regulation.

**Application:** Functional assay

**Figure 1**



**Figure 1.** Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

### References:

- Morfis *et al.* (2008) Receptor Activity-Modifying Proteins Differentially Modulate the G Protein-coupling Efficiency of Amylin Receptors. *Endocrinology* 149(11):5423–5431.
- Hay *et al.* (2005) Pharmacological Discrimination of Calcitonin Receptor: Receptor Activity-Modifying Protein Complexes. *Mol Pharmacol* 67:1655–1665.
- 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.
- RJ Bailey *et al.* (2011). Pharmacological characterization of rat amylin receptors: implications for the identification of amylin receptor subtypes. *BJP* 166:151–167

**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.