

**MULTISCREEN™ DIVISION ARRESTED CELL LINE**  
**HUMAN RECOMBINANT AMY2 (CALCITONIN+RAMP2) RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1510-1a

**Lot Number:** DC1510-1a-112125

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

**Freeze Medium:** Cell Banker 2

**Host cell:** CHO-K1

**Transfection:** Full-length Human CT cDNA (GenBank Accession Number NM\_001742) with FLAG-tag sequence at the N-terminus and Full-length Human receptor activity modifying protein 2 (RAMP2) cDNA (GenBank Accession Number NM\_005854.2) with myc-tag at the C-terminus

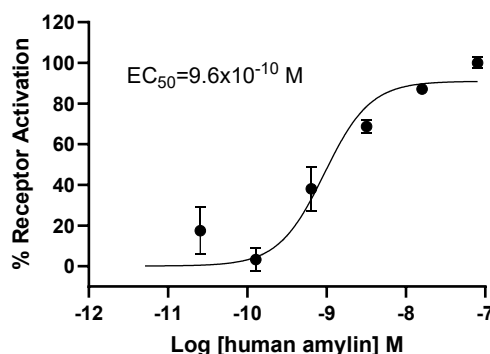
**Recommended Storage:** Liquid nitrogen upon receiving

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

**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 it acts as a high affinity receptor for amylin, a hormone secreted by B cell of pancreas and has a major role in glucose regulation. The Calcitonin/RAMP1 complex has been known as the Amylin subtype AMY1 receptor.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation 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.**

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