

## MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT MAS1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DC1225

**Lot Number:** DC1225-102023

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

**Freeze Medium:** Cell Banker 2  
(Amsbio 11891)

**Host cell:** HEK293T

**Transfection:** Full-length Human MAS1 cDNA (GenBank Accession Number NM\_002377) with FLAG-tag sequence at the N-terminus

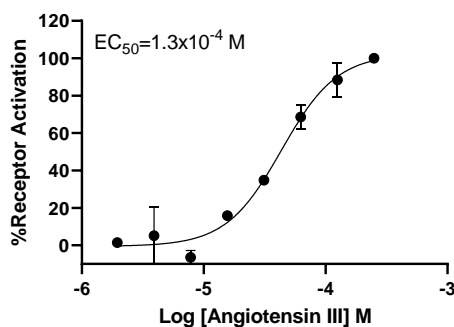
**Recommended Storage:** Liquid nitrogen upon receiving

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

**Background:** MAS1 belongs to the family of MAS-related orphan G protein coupled receptors, which are primarily expressed in primary sensory neurons and mast cells. The MAS-related GPCRs may be implicated in nociception, pruritus, sleep, cell proliferation, circulation, and mast cell degranulation. The MAS1 receptor when activated by binding angiotensin-(1-7) opposes many of the effects of angiotensin-II activated angiotensin receptor. Receptor agonists have similar therapeutic effects as angiotensin-II receptor antagonists including lowering blood pressure. Angiotensin peptide metabolites including angiotensin-III have also been reported to activate MAS1.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

#### References:

Kalyan C. Tirupula<sup>1</sup>, Russell Desnoyer<sup>1</sup>, Robert C. Speth<sup>2</sup>, Sadashiva S. Karnik (2014). Atypical Signaling and Functional Desensitization Response of MAS Receptor to Peptide Ligands. *PLOS ONE*. V9 (7) e103520.

Solinski HJ, Gudermann T, Breit A (2014) Pharmacology and signaling of MAS-related G protein-coupled receptors. *Pharmacol. Rev.* Jul;66(3):570-97.

Young D, Waitches G, Birchmeier C, Fasano O, Wigler M (1986). "Isolation and characterization of a new cellular oncogene encoding a protein with multiple potential transmembrane domains". *Cell*. **45** (5): 711–9.

**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.