

**MULTISCREEN™ STABLE CELL LINE
MOUSE RECOMBINANT AT1 RECEPTOR**

PRODUCT INFORMATION

Catalog Number: Cm1417-1a

Lot Number: Cm1417-1a-012820

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1

Transfection: Expression vector containing full-length mus musculus AGTR1 cDNA (same as GenBank Accession Number NM_177322.3) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DME/F12, 10% FBS, 10 µg/mL puromycin

Stability: Stable for a minimum of 2 months in continuous culture

Data sheet

Background: AT1 (Angiotensin II, type 1) receptor is an angiotensin receptor with vasopressor effects and regulates aldosterone secretion. It is essential in controlling blood pressure and volume in the cardiovascular system. AT1 receptor antagonists are drugs useful in the treatment of hypertension, diabetic nephropathy and congestive heart failure. Blockade of AT1 receptors could also be a novel approach in the treatment of mood disorders and neurodegenerative diseases of the brain.

Application: Functional assays

Figure 1

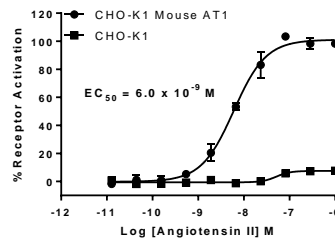


Figure 2

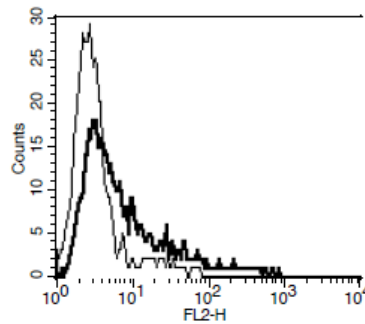


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Bergsma *et al.* (1992) Cloning and characterization of a human angiotensin II type 1 receptor. *Biochem Biophys Res Commun* 183:989-995.

Jagadeesh (1998) Angiotensin II receptors-antagonists, molecular biology, and signal transduction. *Indian J Exp Biol* 36:1171-1194.

Saavedra *et al.* (1998) Blockade of brain angiotensin II AT1 receptors ameliorates stress, anxiety, brain inflammation and ischemia: Therapeutic implications. *Psychoneuroendocrinology* 36:1-18.

FOR RESEARCH USE ONLY.

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