

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT AT1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: H1417

Lot Number: H1417-071211

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 AGTR1 cDNA (same as GenBank Accession Number NM_000685.4 except for one silent mutation) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

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

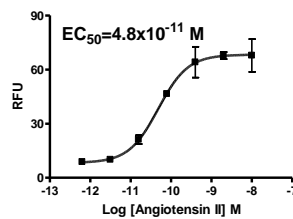


Figure 2

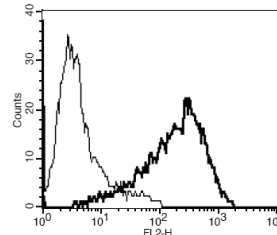


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.