

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT UT RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1035-1a

Lot Number: C1035-1a-090911

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Full-length human UTS2R cDNA (GenBank Accession Number NM_018949.1)

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: In progress

Background: The urotensin 2 receptor (UT or GPR14) is high affinity receptor for urotensin-2 and urotensin-2B. In human, Urotensin II is the most potent vasoconstrictor known, possibly contributing to several human cardiovascular diseases. Both UT receptor and urotensin II are widely expressed in many other tissues, implicating urotensin II in the pathogenesis of a variety of disease processes ranging from hypertension to hepatic cirrhosis and UT antagonists for the possibility of a new range of therapeutic drugs.

Application: Functional assays

Figure 1

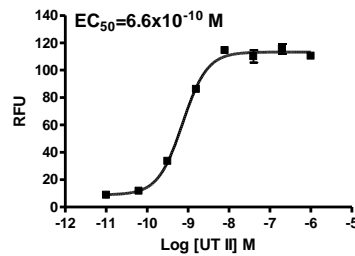


Figure 2

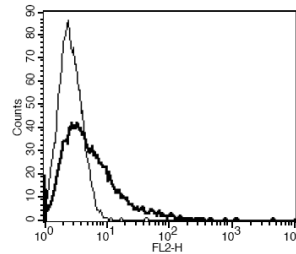


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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:

Kemp *et al.* (2005) Urotensin II: a vascular mediator in health and disease. *Curr Vasc Pharmacol* 3:159-168.

Carotenuto *et al.* (2004) Urotensin-II receptor peptide agonists. *Med Res Rev* 24:577-588.

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.