

## MULTISCREEN™ STABLE CELL LINE RAT RECOMBINANT UT RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1035-1

**Lot Number:** C1035-1-073005

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO dhfr<sup>-</sup>

**Transfection:** Full-length rat Uts2r cDNA (GenBank Accession Number U32673)

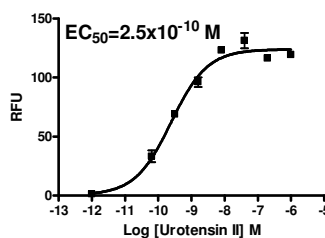
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** Alpha-MEM (Hyclone: SH30265), 10% FBS, 800  $\mu$ g/mL G418

**Stability:** Stable in culture for minimum of two months

**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 legend.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

#### 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.**

© 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.