

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT 5HT4 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1330

**Lot Number:** C1330-012610

**Quantity:** 1 vial (2 x 10<sup>6</sup>) frozen cells

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human HTR4 cDNA (GenBank accession number NM\_000870.4) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

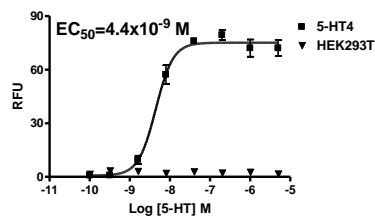
**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin

**Stability:** Stable after minimum two months continuous growth

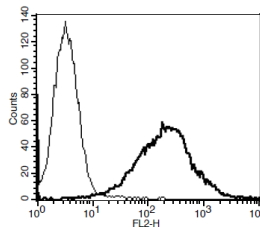
**Background:** The human serotonin receptor 5-HT<sub>4</sub> is a G protein-coupled receptor. 5-HT<sub>4</sub> receptors can be found in the central nervous system and in peripheral tissues such as the alimentary tract, urinary bladder, heart and adrenal gland. 5-HT<sub>4</sub> ligands may prove to be therapeutic in the treatment of disorders such as cardiac arrhythmias, irritable bowel syndrome and urinary incontinence.

**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:

Talley (2003) Evaluation of drug treatment in irritable bowel syndrome. *Br J Clin Pharmacol* 56:362-369.

Van den Wyngaert *et al.* (1997) Cloning and expression of a human serotonin 5-HT<sub>4</sub> receptor DNA. *J Neurochem* 69:1810-1819.

**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.