

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT 5-HT4 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1330-1

Lot Number: 02/20/13

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

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/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: The human serotonin receptor 5-HT4 is a G protein-coupled receptor. 5-HT4 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-HT4 ligands may prove to be therapeutic in the treatment of disorders such as cardiac arrhythmias, irritable bowel syndrome and urinary incontinence.

Application: Functional assays

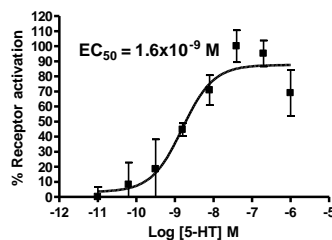


Figure Legend: Dose-dependent stimulation of intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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-HT4 receptor DNA. *J Neurochem* 69:1810-1819.

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.