

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT 5-HT4D RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1520-1

Lot Number: C1520-1-042318

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1

Transfection: Expression vector containing full-length human HTR4 cDNA (GenBank accession number NM_001040172.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 µg/mL puromycin

Stability: Stable in culture for minimum of 1 month

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 1

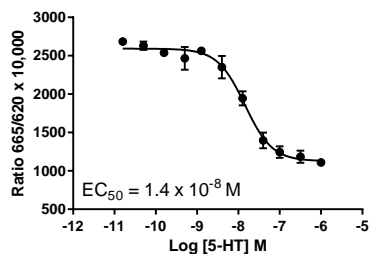


Figure 1. Dose-dependent stimulation of intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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.