

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT 5-HT4B RECEPTOR

PRODUCT INFORMATION

Catalog Number: CG1330

Lot Number: CG1330-022522

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: HEK293T Gqi5

Transfection: Expression vector containing full-length human HTR4 cDNA (GenBank accession number NM_000870.4) with FLAG tag sequence at N-terminus

Propagation Medium: DMEM, 10% FBS, 10 µg/mL puromycin, 50 µg/mL hygromycin

Recommended Storage: Liquid nitrogen upon receiving

Stability: Stable for a minimum of 2 months in continuous culture

Data sheet

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

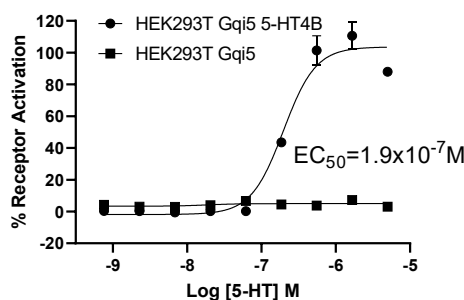


Figure 2

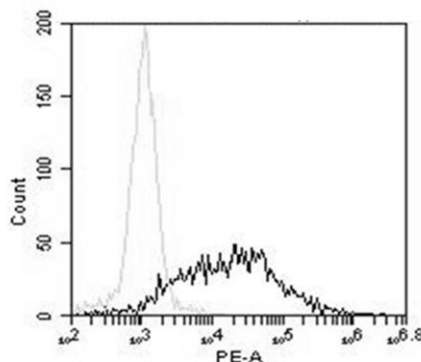


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Black line: parental cells; gray 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-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.