

MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT 5-HT2C RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1326-1

Lot Number: DC1326-1-081222

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: CHO-K1

Transfection: Full-length Human
HTR2C cDNA (GenBank Accession
Number NM_000868) with FLAG-tag
sequence at the N-terminus

Recommended Storage: Liquid
nitrogen upon receiving

Propagation Medium: DMEM/F12,
10% FBS

Background: 5-HT2C (5-hydroxytryptamine receptor 2C) is a receptor for serotonin. It is expressed in the brain and spinal cord, especially the choroid plexus. 5-HT2C receptor agonists may have important clinical value in the treatment of mental and eating disorders, such as depression, panic anxiety, OCD, bulimia and obesity.

Application: Functional assays

Figure 1

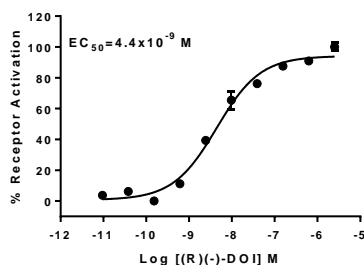


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Martin *et al.* (1998) 5-HT2C receptor agonists: Pharmacological characteristics and therapeutic potential. *J Pharmacol Exp Ther* 286:913-924.

Ni and Miledi (1997) Blockage of 5HT2C serotonin receptors by fluoxetine (Prozac). *Proc Natl Acad Sci USA* 94:2036-2040.

Porter *et al.* (1999) Functional characterization of agonists at recombinant human 5-HT2A, 5-HT2B, and 5-HT2C receptors in CHO-K1 cells. *Br J Pharmacol* 128:13-20.

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.