

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

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1326-1

**Lot Number:** DC1326-1-08/19/09

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

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

**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

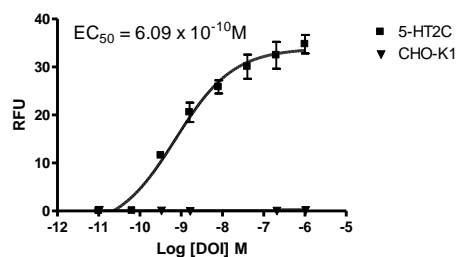
**Propagation Medium:** DMEM/F12, 10% FBS

**Stability:** Stable for 1-2 days after thawing

**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.