

**MULTISCREEN™ STABLE CELL LINE
MOUSE RECOMBINANT 5-HT2A RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: Cm1324-1

Lot Number: Cm1324-1-013023

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1

Transfection: Full-length Mouse HTR2A cDNA (GenBank Accession Number NM_172812) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum two months continuous growth

Background: 5-HT2A (5-hydroxytryptamine receptor 2A) is a receptor for serotonin. It is expressed throughout the central nervous system in the neocortex and olfactory tubercle. Additionally, it is expressed in platelets, fibroblasts and neurons of the peripheral nervous system. 5-HT2A receptor agonists may have important clinical value in the treatment of various disorders, such as depression, anxiety, bipolar disorder and schizophrenia. It is also a receptor for the human polyomavirus, JC virus.

Application: Functional assays

Figure 1

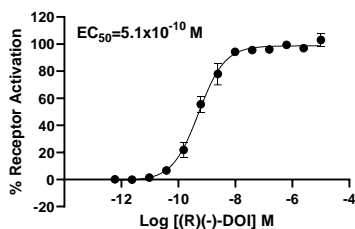


Figure 2

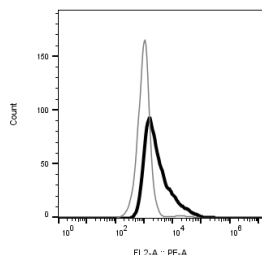


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick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

References:

Meltzer and Li (2003) Serotonin receptors: their key role in drugs to treat schizophrenia. *Prog Neuropsychopharmacol Biol Psychiatry* 27:1159-1172.

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.

Stam *et al.* (1992) Genomic organization, coding sequence and functional expression of human 5-HT2 and HT1A receptor genes. *Eur J Pharmacol* 227:158-162.

FOR RESEARCH USE ONLY.

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.