

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT 5-HT2A RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1324-1

Lot Number: MC1324-1-030222

Quantity: 6.3mg/mL
(10mg, 1mg)

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: CHO-K1

Transfection: Full-length Human HTR2A cDNA (GenBank Accession Number NM_000621) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving; avoid freeze-thaw

Background: 5-HT_{2A} (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-HT_{2A} 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.

Figure 1

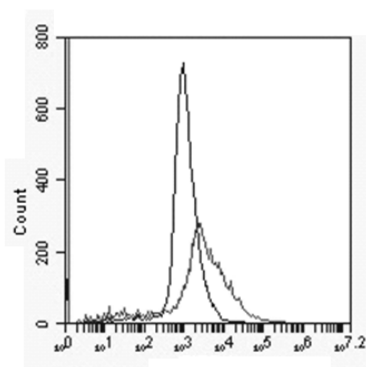


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Black line: parental cells; gray line: receptor-expressing cells.

References:

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-HT_{2A}, 5-HT_{2B}, and 5-HT_{2C} 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 HT_{1A} 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.