

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT 5-HT_{2A} RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1324-1

Lot Number: DC1324-1-053123

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cell Banker 2
(Amsbio)

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

Propagation Medium: DMEM/F12,
10% FBS

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.

Application: Functional assays

Figure 1

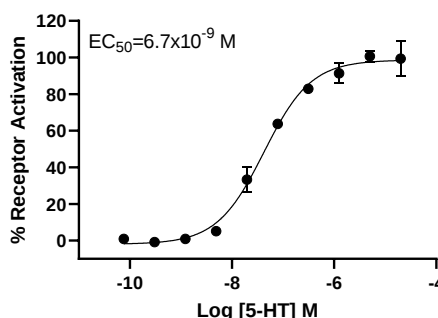


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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.