

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT 5-HT6 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1331-1

Lot Number: DC1331-10/11/2010

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human HTR6 cDNA (GenBank accession number NM_000871) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

Background: Human 5-HT6 receptors are G protein-coupled receptors. The cDNA encodes a 440-amino-acid polypeptide. The 5-HT6 receptors are expressed exclusively in brain regions such as striatum, cortex, nucleus accumbens, olfactory tubercle, hippocampus, hypothalamus, amygdala and cerebellum. The 5-HT6 receptors appear to regulate GABAergic, glutamatergic and cholinergic neurotransmission and are involved in the regulation of cognition, feeding and sleep/wake regulation. They may also play a role in mediating the effects of some antipsychotic agents, and in modulating sensory gating which is often affected in schizophrenic patients.

Application: Functional assays

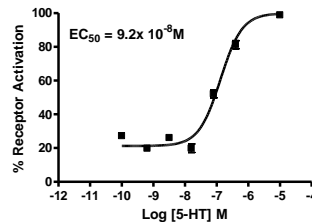


Figure Legend: Dose-dependent increase of intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Wesołowska A (2002) In the search for selective ligands of 5-HT5, 5-HT6 and 5-HT7 serotonin receptors. *Pol J Pharmacol* 54:327-341.

Woolley ML, Marsden CA, and Fone KC (2004) 5-HT6 receptors. *Curr Drug Targets CNS Neurol Disord* 3:59-79.

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.