

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT 5-HT6 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1331-1

Lot Number: C1331-031809

Quantity: 1 vial (2×10^6) 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, 10 μ g/mL puromycin

Stability: Stable after minimum of two months continuous growth

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

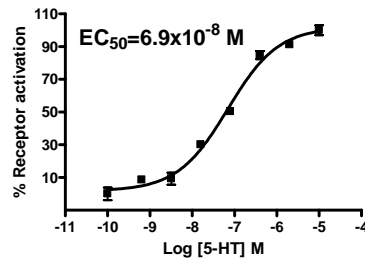


Figure Legend: Dose-dependent increase of intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

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.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.