

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT 5-HT6 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1331-1

**Lot Number:** C1331-031809

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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

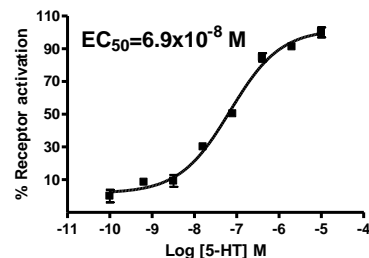
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM/F12, 10% FBS, 10  $\mu$ 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 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.