

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE  
HUMAN RECOMBINANT 5HT7 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1334

**Lot Number:** DC1334-032423

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2  
(Amsbio)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human HTR7 cDNA (GenBank accession number NM\_019860.2) with FLAG tag sequence at N-terminus

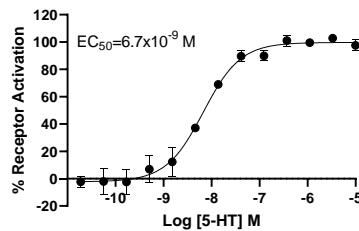
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Background:** The 5-HT7 receptor is a member of the GPCR superfamily of cell surface receptors and is activated by the neurotransmitter serotonin. The 5-HT7 receptor is expressed in a variety of tissues, particularly in the brain, the gastrointestinal tract, and various blood vessels. The 5-HT7 receptor plays a role in smooth muscle relaxation within the vasculature and in the gastrointestinal tract, and is involved in thermoregulation, circadian rhythm, learning and memory, and sleep. This receptor is also involved in mood regulation, suggesting that it may be a useful target in the treatment of depression.

**Application:** Functional assays

**Figure 1**



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

**References:**

Bard *et al.* (1993) Cloning of a novel human serotonin receptor (5-HT7) positively linked to adenylate cyclase. *J Biol Chem* 268:23422-23426.

Lovenberg *et al.* (1993) A novel adenylyl cyclase-activating serotonin receptor (5-HT7) implicated in the regulation of mammalian circadian rhythms. *Neuron* 11:449-458.

**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.