

**MULTISCREEN™ STABLE CELL LINE  
HUMAN RECOMBINANT 5-HT1B RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1320

**Lot Number:** C1320-042110

**Quantity:** 1 vial (2 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human HTR1B cDNA (GenBank accession number NM\_000863.1) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

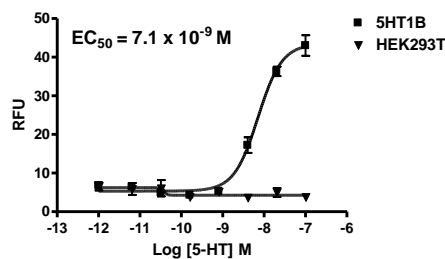
**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin

**Stability:** Stable in culture for minimum of two months

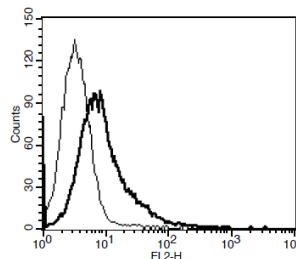
**Background:** The human serotonin receptor 5-HT1B is a G protein-coupled receptor. 5-HT1B receptors are present in many parts of the central nervous system, but most notably, can be found in the basal ganglia, striatum, and frontal cortex. 5-HT1B receptors inhibit the release of many neurotransmitters, such as serotonin, GABA, acetylcholine, and glutamate. 5-HT1B ligands may prove to be therapeutic in the treatment of various disorders such as depression, anxiety, and aggression.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

**References:**

Hamblin *et al.* (1992) Molecular cloning and functional characterization of a human 5-HT1B serotonin receptor: a homologue of the rat 5-HT1B receptor with 5-HT1D-like pharmacological specificity. *Biochem Biophys Res Commun*, 184:752-759.

Hamon *et al.* (1990) The main features of central 5-HT1 receptors. *Neuropsychopharmacology* 3:349-360.

Ruf *et al.* (2009) The 5-HT(1B) receptor: a novel target for the pathophysiology of depression. *Curr Drug Targets* 10:1118-1138.

**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.