

## MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT 5-HT2B RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** Cm1325

**Lot Number:** Cm1325-111513

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

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

**Host cell:** HEK293T

**Transfection:** Full-length Mouse HTR2B cDNA (GenBank Accession Number NM\_008311) with FLAG-tag sequence at the N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

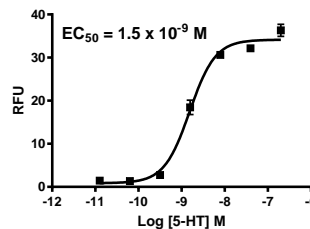
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

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

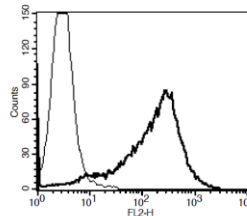
**Background:** 5-HT2B (5-hydroxytryptamine receptor 2B) is a receptor for serotonin. It is expressed in many peripheral and central nervous system tissues, including stomach fundus, liver kidney, muscle, intestine and brain. 5-HT2B receptors are responsible for many cardiovascular and central nervous system functions, such as blood vessel contraction, platelet shape changes, neuronal sensitization to tactile stimuli, and mediation of the hallucinogenic effects of phenylisopropylamin hallucinogens. It has also been shown to be required for heart development.

**Application:** Functional assays

**Figure 1.**



**Figure 2.**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **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:

Bonhaus *et al.* (1995) The pharmacology and distribution of human 5-hydroxytryptamine 2B (5-HT2B) receptor gene products: comparison with 5-HT2A and 5-HT2C receptors. *Br J Pharmacol* 115:622-628.

Nebegil *et al.* (1992) Serotonin 2B receptor is required for heart development. *Proc Natl Acad Sci USA* 97:9508-9513.

Porter *et al.* (1999) Functional characterization of agonists at recombinant human 5-HT2A, 5-HT2B, and 5-HT2C receptors in CHO-K1 cells. *Br J Pharmacol* 128:13-20.

**FOR RESEARCH USE ONLY.**

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