

## MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT LPA5 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** Cm1145-6

**Lot Number:** C1145-6-062012

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

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

**Host cell:** RH7777

**Transfection:** Expression vector containing full-length mouse LPA5 cDNA (GenBank Accession Number: NM\_001163268.1) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

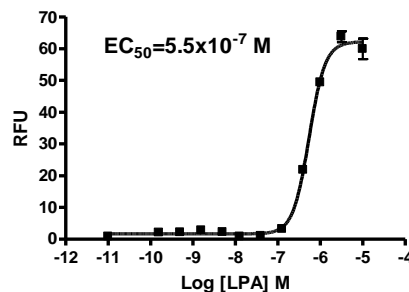
**Propagation Medium:** DMEM, 10% FBS, 3  $\mu$ g/mL puromycin

**Stability:** In progress

**Background:** LPA5 (GPR92) represents a putative orphan G protein coupled receptor with unknown functions. The mouse LPA5 receptor gene has an 81% homology to the human ortholog. Research has shown LPA5 to be expressed in the mouse intestine and upon stimulation with LPA, activates Na<sup>+</sup>/H<sup>+</sup> exchanger 3 and stimulates sodium and fluid absorption. ESTs for LPA5 have been isolated from human normal brain, testis, and tonsil B-cell libraries and from a human blood cancer cell library (pre-B cell). LPA5 protein contains 372 amino acids and shared 36 to 40% sequence identity in the transmembrane regions with the G protein-coupled purinergic receptor P2Y5 (P2RY5), GPR23, and GPR17.

**Application:** Functional assays

**Figure 1**



**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:

Lee *et al.* (2001) Discovery and mapping of ten novel G protein coupled receptor genes. *Gene* 275:83-91.

White *et al.* (2000) Autosomal dominant hypophosphataemic rickets is associated with mutations in FGF23. *Nat Genet* 26:345-348.

Lee *et al.* (2006) GPR92 as a new G12/13- and Gq-coupled lysophosphatidic acid receptor that increases cAMP, LPA5. *J Biol Chem* 281:23589-23597.

Yoo *et al.* (2011) Lysophosphatidic acid 5 receptor induces activation of Na<sup>+</sup>/H<sup>+</sup> exchanger 3 via apical epidermal growth factor receptor in intestinal epithelial cells. *American Journal of Physiology Cell Physiology* Vol. 301 no. 5 C1008-C1016.

**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.