

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT LPA2 (EDG4) RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1050-6

**Lot Number:** C1050-6-061808

**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 human EDG4 cDNA (GenBank Accession Number NM\_004720.4) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

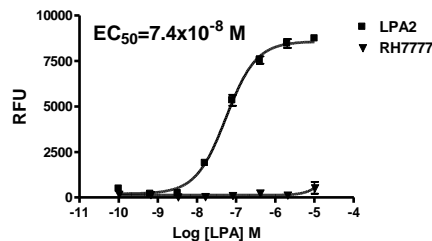
**Propagation Medium:** DMEM, 10% FBS, 5  $\mu$ g/mL puromycin

**Stability:** Stable after minimum of two months continuous growth

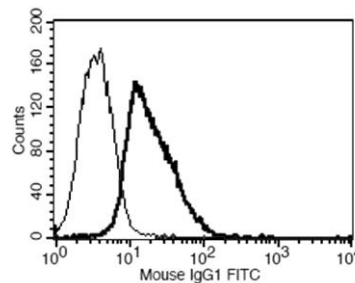
**Background:** The lysophosphatidic acid receptor LPA2 or endothelial differentiation, G-protein coupled receptor 4(EDG-4) is expressed most abundantly in testes and peripheral blood leukocytes. It is reported to be a distinctive functional marker for ovarian carcinoma.

**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:**

Huang *et al.* (2002) Lysophospholipid mediators of immunity and neoplasia. *Biochim Biophys Acta* 1582:161-167.

An *et al.* (1998) Characterization of a novel subtype of human G protein-coupled receptor for lysophosphatidic acid. *J Biol Chem* 273:7906-7910.

**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.