

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

PRODUCT INFORMATION

Catalog Number: C1050-1

Lot Number: C1050-1-082509

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

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-F12, 10% FBS, 10 μ g/mL puromycin

Stability: Stable after minimum of two months continuous growth

Data sheet

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

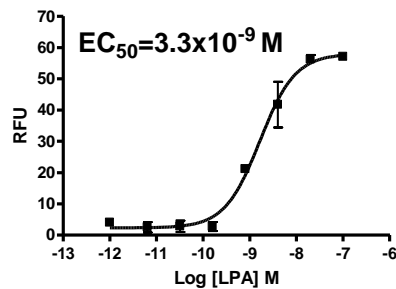


Figure 2

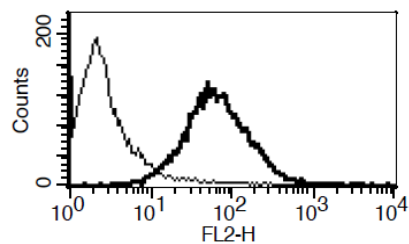


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick 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.