

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

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1050-6

Lot Number: 06/21/13

Quantity: 1 vial (2×10^6) 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

Stability: Stable for 1-2 days after thawing

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

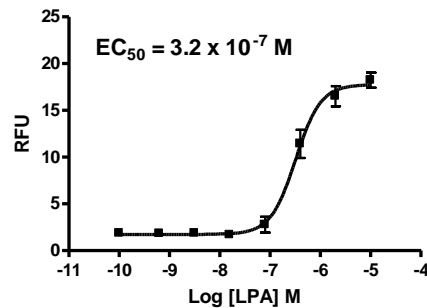


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

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 proteincoupled receptor for lysophosphatidic acid. *J Biol Chem* 273:7906-7910.

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.