

**MULTISCREEN™ MEMBRANE PREPARATION
HUMAN RECOMBINANT LPA3 (EDG7) RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1053-6

Lot Number: MC1053-6-02082012

Quantity: 6.1 mg/ml, 1 mg

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: RH7777

Transfection: Expression vector containing full-length human LPAR3 cDNA (GenBank Accession Number NM_012152) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Background: LPA and the structurally related lysophospholipid mediator sphingosine 1-phosphate (S1P) signal cells through a set of G protein-coupled receptors known as EDG receptors. Some EDG receptors (e.g., EDG1) are S1P receptors; others (e.g., EDG2) are LPA receptors. LPA3 receptor (EDG7) mediates responses preferentially to unsaturated LPA, whereas LPA2 receptor (EDG4) mediates responses to both saturated and unsaturated LPA.

Figure 1

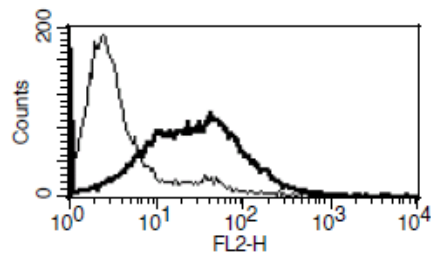


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Bandoh *et al.* (1999) Molecular cloning and characterization of a novel human G-protein-coupled receptor, EDG7, for lysophosphatidic acid. *J. Biol. Chem.* 274:27776-27785.

Ye *et al.* (2005) LPA3-mediated lysophosphatidic acid signalling in embryo implantation and spacing. *Nature* 435:104-108.

FOR RESEARCH USE ONLY.

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