

**MULTISCREEN™**  
**HUMAN ENDOGENOUS LPA1 (EDG-2) RECEPTOR**

---

**PRODUCT INFORMATION**

**Catalog Number:** C1048-E

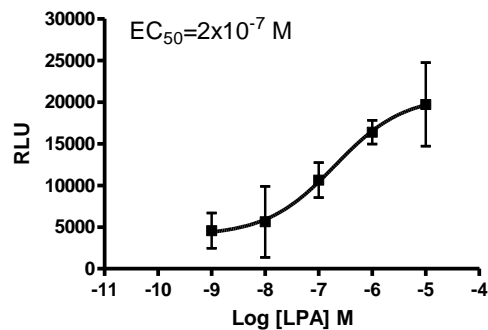
**Lot Number:** C1048-E-083014

**Cell:** Human Fibroblast-Like  
Synoviocytes naturally expressing  
LPA1

**Data sheet**

**Background:** The lipid growth factor lysophosphatidic acid (LPA) is responsible for cell signaling in diverse pathways including survival, proliferation, motility, and differentiation. LPA acts upon target cells by activating its cognate receptors, which belong to the G protein-coupled endothelial differentiation gene (EDG) family. Four mammalian cell surface LPA receptors have been identified so far: EDG-2 (LPA1), EDG-4 (LPA2), EDG-7 (LPA3) and LPA4 (GPR23/P2Y9).

**Application:** Functional assays



**Figure legend:** Dose-dependent stimulation of cell migration upon treatment with ligand.

**References:**

Mills and Moolenaar (2003) The emerging role of lysophosphatidic acid in cancer. *Nat Rev Cancer* 3:582-591.

Yang *et al.* (2002) In vivo roles of lysophospholipid receptors revealed by gene targeting studies in mice. *Biochim Biophys Acta* 1582:197-203.

**FOR SCREENING BY MULTISPAN INC. ONLY.**  
**NOT SOLD AS PRODUCT.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

26219 Eden Landing Road, Hayward, California 94545-3718 Phone: (510) 887-0817 info@multispaninc.com