

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT LPA4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1087-1a

Lot Number: C1087-1a-040811

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 LPA4 cDNA (GenBank accession number NM_005296) 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 one month continuous growth

Background: Human LPA4 (also known as P2Y9 and GPR23) receptors are G protein-coupled receptors. The cDNA encodes a 370-amino-acid polypeptide. The LPA4 receptors are predominately expressed in ovary, and also present in kidney, skeletal muscle and other brain and peripheral tissues. In rat neuroblastoma cells overexpressing the receptors, LPA promotes neurite retraction and cell rounding indicating LPA4 may mediate morphological changes in neuronal cells. Studies in LPA4 knockout mice revealed a role of LPA4 in the negative control of cell migration.

Application: Functional assays

Figure 1

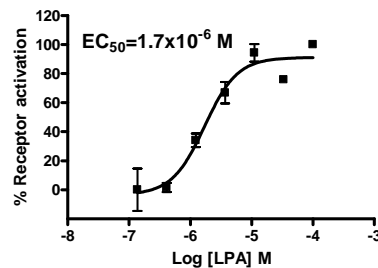


Figure 2

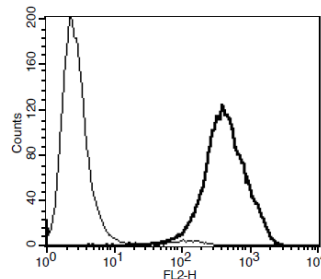


Figure 1. Dose-dependent increase of intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

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:

Yanagida *et al* (2007) LPA4/p2y9/GPR23 mediates rho-dependent morphological changes in a rat neuronal cell line. *J Biol Chem* 282:5814-5824.

Lee *et al* (2008) Role of LPA4/p2y9/GPR23 in negative regulation of cell motility. *Mol Biol Cell* 19:5435-5445.

Noguchi *et al.* (2003) Identification of p2y9/GPR23 as a novel G protein-coupled receptor for lysophosphatidic acid, structurally distant from the EDG family. *J Biol Chem* 278:25600-25606.

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.