

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT GPR109B RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1059-1

Lot Number: C1059-1-040709

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 GPR109B cDNA (GenBank Accession Number NM_006018.1) 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 two months of continuous growth

Background: GPR109B (or HM74) is a 387-amino acid, 7-transmembrane protein responsible for the directed migration of specific cell types. GPR109B has highest expression in spleen, lymphocytes and adipose tissue, and lower expression in the heart, placenta, prostate and bone marrow. Nicotinic acid has been used as a lipid-lowering agent possibly by acting through GPR109A and GPR109B in adipose tissues. This supports the suggestion that signaling through this nicotinic acid receptors may be a mechanism for the treatment of hyperlipidemia and insulin-resistant states.

Application: Functional assays

Figure 1

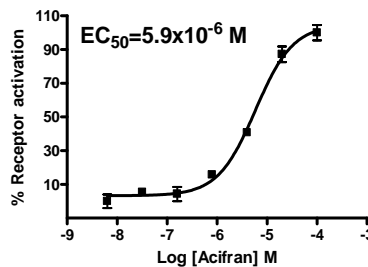


Figure 2

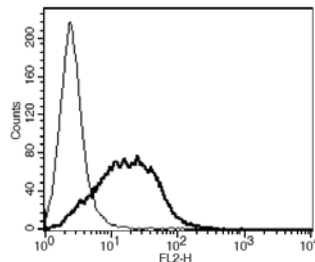


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Karpe and Frayn (2004) The nicotinic acid receptor-a new mechanism for an old drug. *Lancet* 363:1892-1894.

Tunaru *et al.* (2003) PUMA-G and HM74 are receptors for nicotinic acid and mediate its anti-lipolytic effect. *Nature Med* 9:352-355.

Wise *et al.* (2003) Molecular identification of high and low affinity receptors for nicotinic acid. *J Biol Chem* 278:9869-9874.

Nomura *et al.* (1993) Molecular cloning of cDNAs encoding a LD78 receptor and putative leukocyte chemotactic peptide receptors. *Int Immun* 5:1239-1249.

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.