

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR119 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1298

Lot Number: C1298-081211

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human GPR119 cDNA (GenBank Accession Number NM_178471.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: In progress

Background: GPR119 is a G protein-coupled receptor for the endogenous lipid signaling agent oleoylethanolamide (OEA). GPR119 is predominantly expressed in the pancreas (beta-cells) and gastrointestinal tract (enteroendocrine cells). In vitro studies have indicated a role for GPR119 in the modulation of insulin release by pancreatic beta-cells and of GLP-1 secretion by gut enteroendocrine cells.

Application: Functional assays

Figure 1

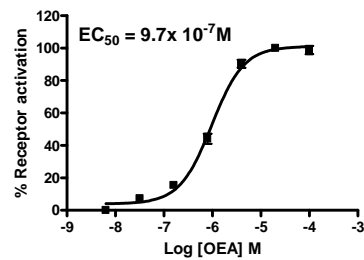


Figure 2

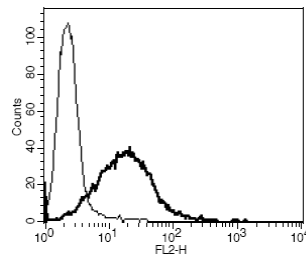


Figure Legend: **Figure 1.** Dose-dependent accumulation of intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells;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

References:

Hilary *et al.* (2006) Deorphanization of a G protein-coupled receptor for oleoylethanolamide and its use in the discovery of smallmolecule hypophagic agents. *Cell Metabolism* 3:167-175.

Overton *et al.* (2007) GPR 119, a novel G protein-coupled receptor target for the treatment of type 2 diabetes and obesity. *Br J Pharmacol* 153 suppl 1:S76-S81.

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.