

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR119 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1298-1a

**Lot Number:** C1298-1a-121009

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human GPR119 cDNA (GenBank Accession Number NM\_178471.1) with FLAG tag sequence at N-terminus

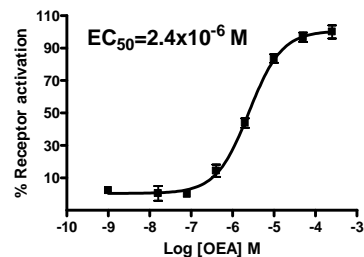
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM-F12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** Stable in culture for minimum of two months

**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 Legend:** Dose-dependent accumulation of intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

#### 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.