



**MULTISPAN**  
Assay Experts in GPCRs and Beyond

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT GPR43 RECEPTOR**

**PRODUCT INFORMATION**

**Catalog Number:** DC1104

**Lot Number:** DC1104-020924

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker2

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human FFAR2 cDNA (GenBank Accession Number NM\_005306) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

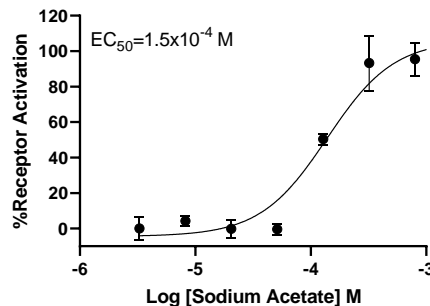
**Propagation Medium:** DMEM, 10% FBS

**Data Sheet**

**Background:** GPR43, or free fatty acid receptor 2 (FFAR2), encodes a deduced 330-amino acid protein with 7 transmembrane domains. GPR43 is expressed by enteroendocrine L cells containing peptide YY in the human large intestine. The receptor binding of short-chain fatty acids potentially provides a molecular link between diet, gastrointestinal bacterial metabolism, and immune and inflammatory responses.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

**References:**

Sawzdargo et al. (1997) Cluster of four novel human G protein-coupled receptor genes occurring in close proximity to CD22 gene on chromosome 19q13.1. *Biochem Biophys Res Commun* 239:543-547.

**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.

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