

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

**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. The receptor binding of short-chain fatty acids potentially provides a intestine molecular link between diet, gastrointestinal bacterial metabolism, and immune and inflammatory responses.

**PRODUCT INFORMATION**

**Catalog Number:** DC1104

**Lot Number:** DC1104-012412

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

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

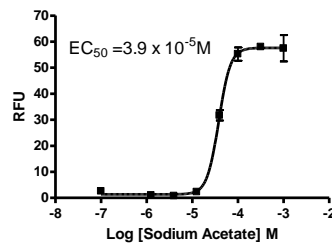
**Host cell:** HEK293T

**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

**Stability:** stable 1-2 days after thawing



**Figure 1.** Dose-dependent 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.

**Application:** Functional assays

**Figure 1**

**FOR RESEARCH USE ONLY.**

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.