

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
MOUSE RECOMBINANT GPBAR1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DCm1361-1

Lot Number: 06/28/11

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Expression vector containing full-length mouse GPBAR1 cDNA (GenBank Accession Number BC116914) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS,

Stability: Stable for 1-2 days after thawing

Background: The G protein-coupled bile acid receptor GPBAR1 (or GPR131) mediates bile acids-induced rapid elevation of intracellular cAMP levels. It is implicated in the suppression of macrophage functions and regulation of energy homeostasis by bile acids.

Application: Functional assays

Figure 1

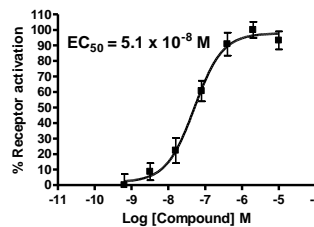


Figure 1. Dose-dependent increase of intracellular cAMP upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Maruyama *et al.* (2002) Identification of membrane-type receptor for bile acids (M-BAR). *Biochem Biophys Res Commun* 298:714-719.

Katsuma *et al.* (2005) Bile acids promote glucagon-like peptide-1 secretion through TGR5 in a murine enteroendocrine cell line STC-1. *Biochem Biophys Res Commun* 329:386-390.

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.