

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT GPR84 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DCG1137-1

Lot Number: DCG1137-1-112625

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1 Gαq15

Transfection: Expression vector containing full-length human GPR84 cDNA (GenBank accession Number NM_020370) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DME/F12, 10% FBS

Background: Human GPR84 is a G protein-coupled receptor with 396 amino acids. The hGPR84 transcript was found in brain, heart, muscle, colon, thymus, spleen, kidney, liver, intestine, placenta, lung and leukocytes. A study in GPR84-deficient mice revealed a novel role for GPR84 in regulating early IL-4 gene expression in activated T cells in response to CD3 crosslinking.

Application: Functional assays

Figure 1

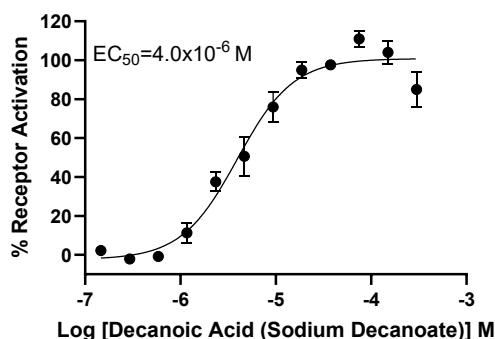


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with MULTISCREEN™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Wittenberger *et al.* (2001) An expressed sequence tag (EST) data mining strategy succeeding in the discovery of new G-protein coupled receptors. *J Mol Biol* 307:799-813.

Venkataraman and Kuo (2005) The G-protein coupled receptor, GPR84 regulates IL-4 production by T lymphocytes in response to CD3 crosslinking. *Immunol Lett* 101:144-153.

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.