

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR34 RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** CG1095

**Lot Number:** CG1095-022718

**Quantity:** 1 vial (2 x 10<sup>6</sup>) frozen cells

**Freeze Medium:** Cellbanker 2 (Amsbio 11891)

**Host cell:** HEK293T Gqi5

**Transfection:** Expression vector containing full-length Human GPR34 cDNA (GenBank Accession Number: NM\_005300) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin, 250 µg/mL hygromycin

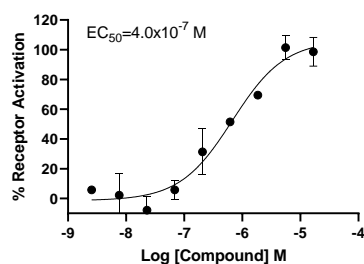
**Stability:** In progress

### Data sheet

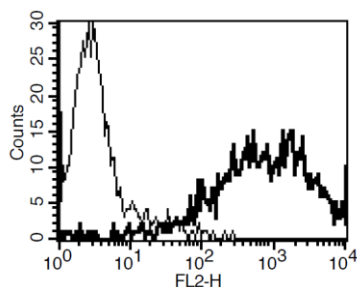
**Background:** GPR34 is an orphan G protein coupled receptor, belongs to P2Y receptor family. It is a potent mediator of lysophospholipid and expressed in various tissues including stomach Mast cells, placenta, spleen, thymus, ovary, Heart, brain etc. GPR34 receptor is responsible for enhancing the antigen induced degranulation of mast cells.

**Application:** Functional assays

**Figure 1**



**Figure 2**



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

**Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

### References:

- K. Makide and J. Aoki (2013) GPR34 as a lysophosphatidylserine receptor. *J. Biochem.* 2013;153 (4):327–329.
- H. Kitamura *et al.* (2012) GPR34 is a receptor for lysophosphatidylserine with a fatty acid at the sn-2 position *J. Biochem.* 2012;151 (5):511–518
- T. Sugo *et al.* (2006) Identification of a lysophosphatidylserine receptor on mast cells. *Biochemical and Biophysical Research Communications* 341 (2006) 1078–1087

**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.