

**MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT GPR34 RECEPTOR**

PRODUCT INFORMATION

Catalog Number: DCG1095

Lot Number: DCG1095-050818

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

Freeze Medium: Cellbanker 2 (Amsbio 11891)

Host cell: HEK293T Gαq_{i5}

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

Stability: 1-2 days after thawing

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

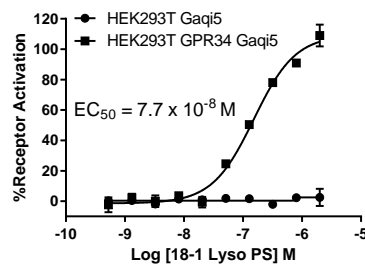


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.