

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR15 RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1069

Lot Number: C1069-012722

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: HEK293T

Transfection: Expression vector containing full-length human GPR15 cDNA (GenBank Accession Number NM_005290.4)

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

Recommended Storage: Liquid nitrogen upon receiving

Stability: Stable in continuous culture for a minimum of 2 months

Data sheet

Background: GPR15 is a class A orphan G protein-coupled receptor found in epithelial cells, synovial macrophages, endothelial cells and lymphocytes especially T cells. Studies have accumulated evidence that GPR15 may modulate chronic inflammation. Findings also have implicated GPR15 for potential therapeutic intervention.

Application: Functional assay

Figure 1

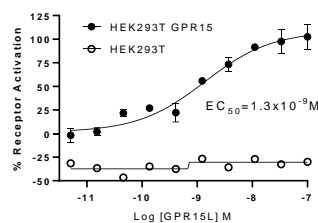


Figure 2

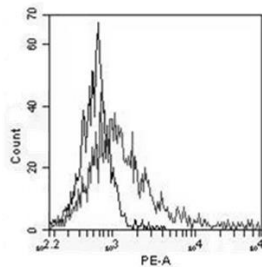


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Ocón, B., Pan, J., Dinh, T. T., Chen, W., Ballet, R., Bscheider, M., Habtezion, A., Tu, H., Zabel, B. A., & Butcher, E. C. (2017). A mucosal and cutaneous chemokine ligand for the lymphocyte chemoattractant receptor GPR15. *Frontiers in Immunology*, 8.

Suply, T., Hannedouche, S., Carte, N., Li, J., Grosshans, B., Schaefer, M., Raad, L., Beck, V., Vidal, S., Hiou-Feige, A., Beluch, N., Barbieri, S., Wirsching, J., Lageyre, N., Hillger, F., Debon, C., Dawson, J., Smith, P., Lannoy, V., Bassilana, F. (2017). A natural ligand for the orphan receptor GPR15 modulates lymphocyte recruitment to epithelia. *Science Signaling*, 10(496).

Hayn, M., Blötz, A., Rodríguez, A., Vidal, S., Preising, N., Ständker, L., Wiese, S., Stürzel, C. M., Harms, M., Gross, R., Jung, C., Kiene, M., Jacob, T., Pöhlmann, S., Forssmann, W.-G., Münch, J., Sparrer, K. M., Seuwen, K., Hahn, B. H., & Kirchhoff, F. (2021). Natural cystatin C fragments inhibit GPR15-mediated HIV and SIV infection without interfering with GPR15L signaling. *Proceedings of the National Academy of Sciences*, 118(3).

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.