

MULTISCREEN™ β -ARRESTIN STABLE CELL LINE HUMAN RECOMBINANT GPR132 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CA1066

Lot Number: CA1066-081619

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Cellbanker 2

Host cell: HEK293T

Transfection: Expression vector containing full-length human GPR132 cDNA (GenBank Accession Number AF083955) with FLAG tag sequence at N-terminus and ARRB2 cDNA (GenBank Accession Number NM_004313.3)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin, 250 μ g/mL hygromycin

Stability: In progress

Background: GPR132 (also known as G2A) is a proton-sensing receptor that is broadly expressed in the appendix, lymph nodes and other tissues. It may also be a receptor for oxidized free fatty acids. It appears to play a role in regulating certain cells of the autoimmune system by responding to the extracellular acidification characteristic of tissues suffering from malignant cell growth or inflammation

Application: Functional assays

Figure 1

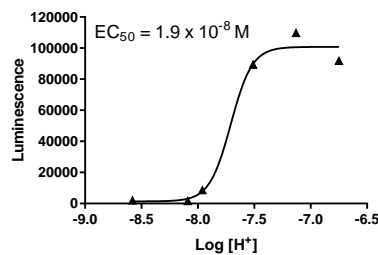


Figure 2

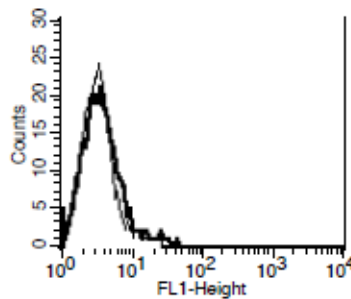


Figure 1. Dose-dependent stimulation from arrestin recruitment upon treatment with ligand, measure with MultiScreen™ β -Arrestin Assay Kit (Multispan MSBAK01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Obinata, *et al.* (2005) Identification of 9-Hydroxyoctadecadienoic Acid and Other Oxidized Free Fatty Acids as Ligands of the G Protein-coupled Receptor G2A. *JOURNAL OF BIOLOGICAL CHEMISTRY* VOL. 280, NO. 49, pp. 40676–40683,

Murakami *et al.* (2004) G2A Is a Proton-sensing G-protein-coupled Receptor Antagonized by Lysophosphatidylcholine. *THE JOURNAL OF BIOLOGICAL CHEMISTRY* Vol. 279, No. 41, Issue of October 8, pp. 42484–42491,

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.