

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR142 RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** C1286a

**Lot Number:** C1286a-120117

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

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human GPR142 cDNA (GenBank Accession Number NM\_181790.1) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

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

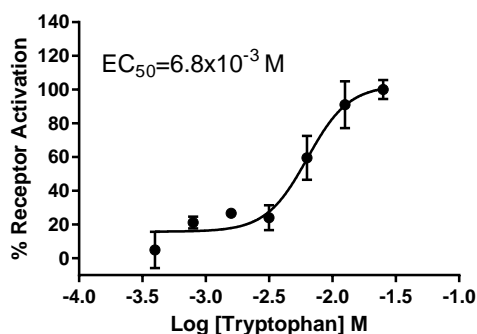
**Stability:** Stable for a minimum of 2 months in continuous culture

### Data sheet

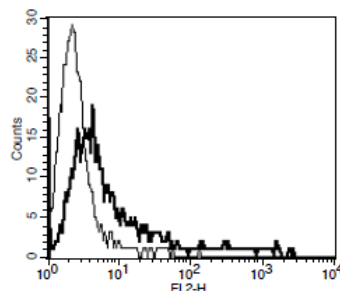
**Background:** GPR142 belongs to the family of class A (rhodopsin-like) orphan G protein-coupled receptors. Research has linked GPR142 to type 2 diabetes mellitus. GPR142 is reported to be highly expressed in pancreatic β-cells, and upon ligand binding and also in the presence of a high concentration of blood glucose, can stimulate insulin secretion. It has been hypothesized that GPR142 agonists can provide a benefit over existing type 2 diabetes therapies because of a greatly reduced risk of hypoglycemia.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose response of intracellular IP1 accumulation upon treatment with ligand, measured with IP-one Tb kit. **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:

Fredriksson R et al. (2003). "Seven evolutionarily conserved human rhodopsin G protein coupled receptors lacking close relatives". *FEBS Lett* 554(3):381-8.

Lizarzaburu M et al. (2012). "Discovery and optimization of a novel series of GPR142 agonists for the treatment of type 2 diabetes mellitus". *Bioorganic & Medicinal Chemistry Letters*

**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.