

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT PAF RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1218

Lot Number: C1218-091505

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human PTAFR cDNA (GenBank Accession Number NM_000952) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum of two months continuous growth

Data sheet

Background: The platelet-activating factor (PAF) receptor mediates a wide range of biological responses to PAF, a potent glycerophospholipid released from a variety of cell types such as stimulated basophils, platelets, polymorphonuclear neutrophils and macrophages. PAF is involved in a diverse array of biological activities related to inflammatory and immune responses as well as cardiovascular, respiratory and nervous system physiology. In humans, various diseases have been associated with PAF, such as allergic asthma, endotoxic shock, atherosclerosis and psoriasis.

Application: Functional assays

Figure 1

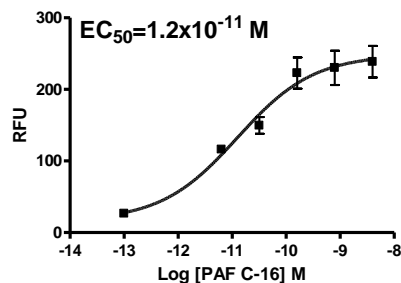


Figure 2

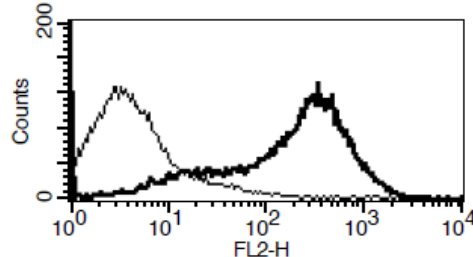


Figure 1. Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Dupre *et al.* (2003) Trafficking, ubiquitination, and down-regulation of the human platelet-activating factor receptor. *J Biol Chem* 278:48228-48235.

Seyfried *et al.* (1992) The human platelet-activating factor receptor gene (PTAFR) contains no introns and maps to chromosome 1. *Genomics* 13:832-834.

Van Biesen *et al.* (1996) G(o)-protein alpha-subunits activate mitogenactivated protein kinase via a novel protein kinase C-dependent mechanism. *J Biol Chem* 271:1266-1269.

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.