

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT P2Y₁ RECEPTOR

PRODUCT INFORMATION

Catalog Number: MC1160-3

Lot Number: MC1160-3-011426

Quantity: 5 mg per vial

Protein concentration: 3.6 mg/mL

Packaging Buffer: 20mM Gly-Gly, 1 mM MgCl₂, 25mM Sucrose (pH 7.2)

Host cell: 1321N1

Transfection: Full-length human P2Y₁ cDNA (GenBank Accession Number NM_002563)

Recommended Storage: Liquid nitrogen upon receiving

Data sheet

Background: P2Y₁ is a receptor for ATP and ADP. P2Y₁ contributes to platelet shape change. A number of P2Y₁ receptor-specific antagonists, such as, MRS-2179, A3P5P, A3P5P and A2P5P have been discovered. They inhibit calcium ion mobilization and shape change in platelets. P2Y₁-deficient mice and mice treated with the P2Y₁ antagonist MRS2179 displayed significantly less arterial thrombosis than their respective controls. Combination of P2Y₁ deficiency with P2Y₁₂ inhibition led to a significant additive effect.

Application: Functional assays

Figure 1

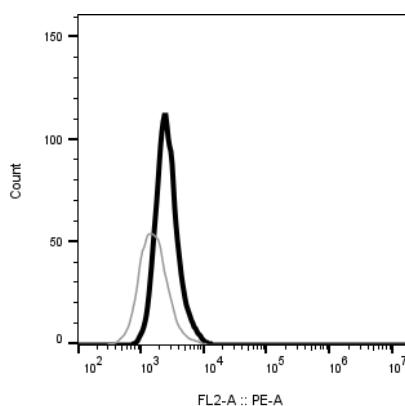


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Baurand and Gachet (2003) The P2Y₁ receptor as a target for new antithrombotic drugs: a review of the P2Y₁ antagonist MRS-2179. *Cardiovasc Drug Rev* 21:67-76.

Jin *et al.* (1998) Molecular basis for ADP-induced platelet activation. II. The P2Y₁ receptor mediates ADP-induced intracellular calcium mobilization and shape change in platelets. *J Biol Chem* 273:2030-2034.

Lenain *et al.* (2003) Inhibition of localized thrombosis in P2Y₁-deficient mice and rodents treated with MRS2179, a P2Y₁ receptor antagonist. *J Thromb Haemost* 1:1144-1149.

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.

26219 Eden Landing Road, Hayward, California 94545-3718 Phone: (510) 887-0817 info@multispaninc.com