

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT P₂Y₁ RECEPTOR**

PRODUCT INFORMATION

Catalog Number: C1160-3

Lot Number: C1160-3-022609

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: 1321N1

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

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 100 µg/mL G418

Stability: Stable after minimum of two months continuous growth

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, A3P5PS, 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

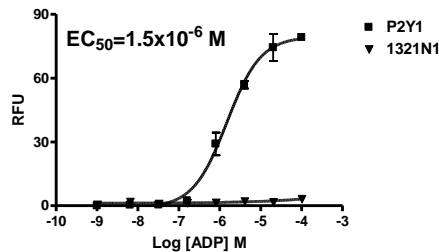


Figure legend: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Baurand and Gachet (2003) The P2Y(1) receptor as a target for new antithrombotic drugs: a review of the P2Y(1) 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.