

MULTISCREEN™ STABLE CELL LINE
RAT RECOMBINANT P2Y12 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1170-1

Lot Number: C1170-1-102008

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO dhfr⁻

Transfection: Full length rat R2RY12 cDNA (GenBank Accession Number N_c_022800)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: α-MEM, 10% FBS, 100 µg/ml hygromycin

Stability: Stable after minimum of two months continuous growth

Background: P2RY12 or P2Y12 is a G-protein coupled receptor for ADP and ATP. ADT released from platelets acts in a positive feedback mechanism on P2Y12 and P2Y1 on platelets and is critical for sustained aggregation and stabilization of thrombin. P2Y12 receptor plays a central role in platelet activation and is the target of antithrombotic drugs such as ticlopidine and clopidogrel.

Application: Functional assays

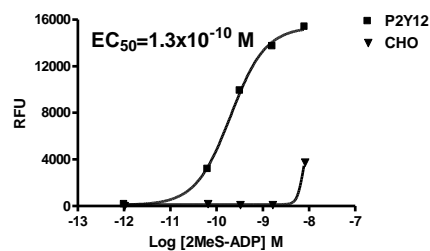


Figure legend: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Hardy *et al.* (2004) Reciprocal cross-talk between P2Y1 and P2Y12 receptors at the level of calcium signaling in human platelets. *Blood* 104:1745-1752.

Cattaneo *et al.* (2003) Molecular bases of defective signal transduction in the platelet P2Y12 receptor of a patient with congenital bleeding. *Proc Natl Acad Sci USA* 100: 1978-1983.

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.