

## MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT P2Y11 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DC1169-3

**Lot Number:** DC1169-3-043021

**Quantity:** 1 vial (4 x 10<sup>6</sup>) frozen cells

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

**Host cell:** 1321N1

**Transfection:** Expression vector containing full-length human P2RY11 cDNA (GenBank Accession Number NM\_002566) with FLAG tag sequence at N-terminus

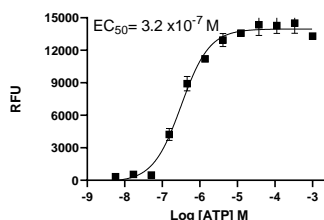
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Background:** P2Y purinoceptor 11 encodes P2Y11 (P2Y11) which is a receptor for ATP and ADP coupled to G-proteins that activate both phosphatidylinositol-calcium and adenylyl cyclase second messenger systems. P2Y11 receptor is unique amongst P2Y receptors with its coupling to the adenylyl cyclase pathway. Through this mechanism, released nucleotides can increase in P2Y11-containing cells. P2Y11 receptor has been reported to be involved in the granulocytic differentiation of promyelocytes and in the maturation of monocyte-derived dendritic cells.

**Application:** Functional assays

**Figure 1**



**Figure 1:** Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01)

#### References:

Torres *et al.* (2002) P2Y11 receptors activate adenylyl cyclase and contribute to nucleotide-promoted cAMP formation in MDCK-D(1) cells. A mechanism for nucleotide-mediated autocrine-paracrine regulation. *J Biol Chem* 277:7761-7765.

Schnurr *et al.* (2003) ATP gradients inhibit the migratory capacity of specific human dendritic cell types: implications for P2Y11 receptor signaling. *Blood* 102:613-620.

**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.