

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT P2Y6 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1164-3

**Lot Number:** C1164-3-040611

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** 1321N1

**Transfection:** Expression vector containing full-length human P2Y6 cDNA (GenBank Accession Number NM\_004154) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

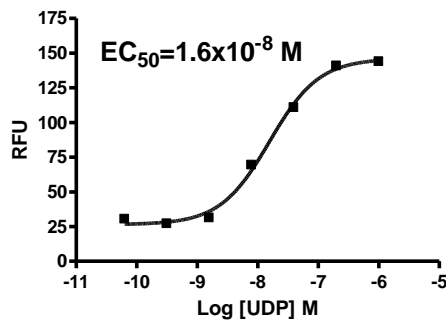
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stable after minimum of two months continuous growth

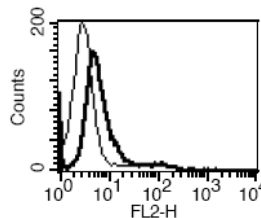
**Background:** The human P2Y6 receptor is a member of the G protein-coupled P2 receptor family that responds specifically to the extracellular nucleotide uridine diphosphate (UDP). The distribution of the P2Y6 receptor is widespread including heart, blood vessels and brain. P2Y6 receptors are significantly upregulated in T cells infiltrating regions of inflamed bowel and are involved in monocytic release of IL-8 in response to UDP or lipopolysaccharide (LPS). The immune effect of UDP/P2Y6 signaling on the production of proinflammatory cytokines has been shown to be selective and dependent on cell types.

**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:

Communi *et al.* (1996) Cloning, functional expression and tissue distribution of the human P2Y6 receptor. *Biochem Biophys Res Commun* 222:303-308.

Warny *et al.* (2001) P2Y(6) nucleotide receptor mediates monocyte interleukin-8 production in response to UDP or lipopolysaccharide. *J Biol Chem* 276:26051-26056.

**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.