

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT A2B RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1429

Lot Number: C1429-080610

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human ADORA2B cDNA (GenBank accession number NM_000676.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: A2B is a receptor for adenosine. A2B receptor is upregulated during intestinal inflammation and mediates key events such as chloride, IL-6 and fibronectin secretion in intestinal epithelial cells. A2B receptor antagonists may have important clinical value in the treatment of inflammatory diseases, such as asthma and chronic obstructive pulmonary disease (COPD), as well as inflammatory bowel disease.

Application: Functional assays

Figure 1

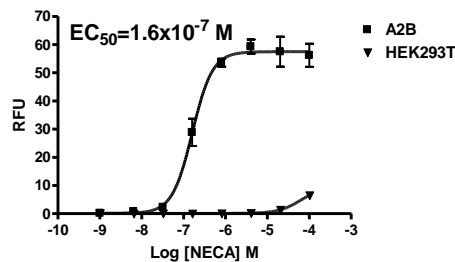


Figure 2

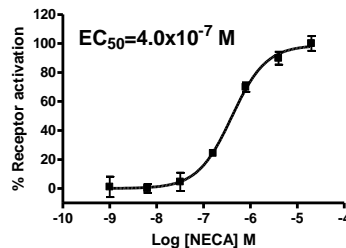


Figure 3

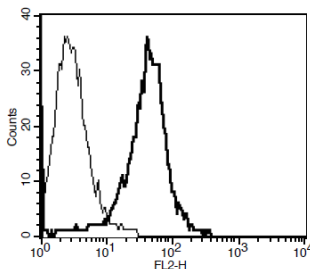


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

Figure 2. Dose-dependent stimulation of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

Figure 3. 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:

Hasko *et al.* (2008) Adenosine receptors: therapeutic aspects for inflammatory and immune diseases. *Nat Rev Drug Discov* 7:759-770.

Linden *et al.* (1999) Characterization of human A (2B) adenosine receptors: radioligand binding, wester blotting, and coupling to G(q) in human embryonic kidney 293 cells and HMC-1 mast cells. *Mol Pharmacol* 56:705-713.

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.