

**MULTISCREEN™ MEMBRANE PREPARATION
RAT RECOMBINANT A2B RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: MCr1429

Lot Number: MCr1429-100818

Quantity: 1 vial (15.7mg/ml, 5mg)

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 250 mM Sucrose

Host cell: HEK293T

Transfection: Expression vector containing full-length rat ADORA2B cDNA (GenBank accession number NM_017161.1) with FLAG tag sequence at N-terminus

Recommended Storage: -80°C upon receiving, avoid repeated freeze-thaw

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. In rats, the A2B receptor has also been shown to stimulate the growth of arterial endothelial cells and may be critical in regulating vascular remodeling associated with endothelial cell proliferation in angiogenesis, collateral vessel development, and recovery after vascular injury.

Application: Functional assays

Figure 1

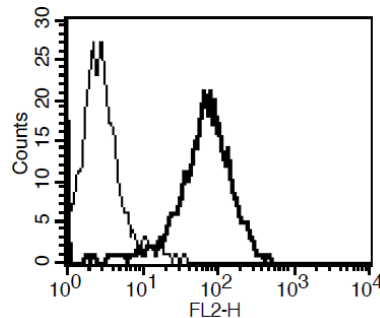


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.

Dubey *et al.* (2002) A2B Adenosine Receptors Stimulate Growth of Porcine and Rat Arterial Endothelial Cells. *Hypertension* 39: 530-535.

FOR RESEARCH USE ONLY.

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.