

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT A3 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1430

Lot Number: C1430-080510

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length Human ADORA3 cDNA (GenBank Accession NM_000677.2) with FLAG-tag sequence at the 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: Adenosine A3 receptors are involved in a variety of intracellular signaling pathways and physiological functions. They are expressed in a wide range of human tissues, but most predominantly in the lung and liver. Recent animal model studies have shown that A3 receptors play important roles in brain ischemia, immunosuppression, and bronchospasm. A3 receptor agonists and/or agonists may have important clinical value in the treatment of asthma and inflammation.

Application: Functional assays

Figure 1

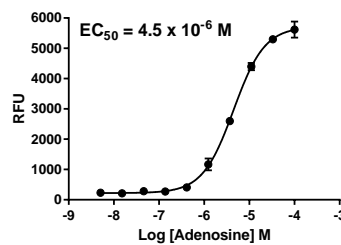


Figure 2

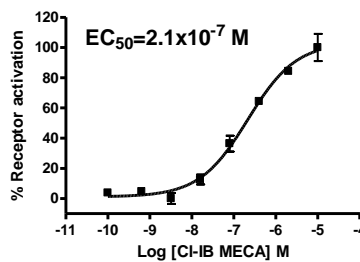


Figure 3

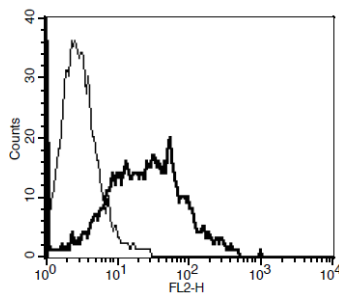


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ 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:

Baraldi *et al.* (2000) A (3) adenosine receptor ligands: history and perspectives. *Med Res Rev* 20:103-128.

Salvatore *et al.* (1993) Molecular cloning and characterization of the human A3 adenosine receptor. *Proc Natl Acad Sci USA* 90:10365-10369.

Schepp *et al.* (2008) Bench-to bedside review: adenosine receptors - promising targets in acute lung injury? *Crit Care* 12:226.

FOR RESEARCH USE ONLY.

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