

**MULTISCREEN™ MEMBRANE PREPARATION
HUMAN RECOMBINANT α2A RECEPTOR**

PRODUCT INFORMATION

Catalog Number: MC1434-1a

Lot Number: MC1434-1a-06192012

Quantity: 1 vial (15.6 mg/ml)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human α2a cDNA (GenBank Accession Number: NM_000681) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen

Data sheet

Background: Adrenergic α2a receptor belongs to the group of nine adrenoceptors that mediate the biological actions of adrenaline and noradrenaline. The α2a receptor is expressed in the brain, spleen, kidney, heart and gastrointestinal tract. It is centrally predominant and exerts a tonic sympathoinhibitory function, whereas peripherally it has a vasoconstrictive effect. The α2a receptor is also involved in sedation, digestive functions and analgesia.

Application: [³⁵S]GTPγS Binding Assay

Figure 1

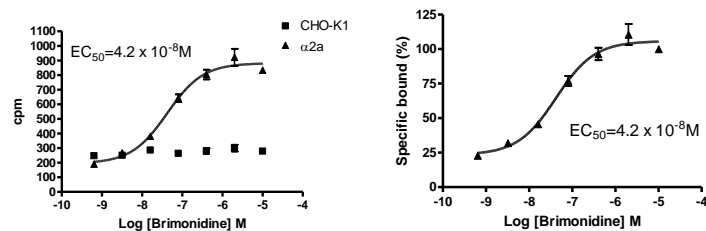


Figure 2

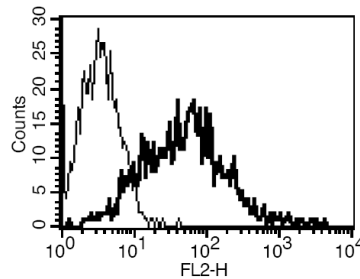


Figure 1. Dose-dependent [³⁵S]GTPγS Binding to α2a membrane upon treatment with agonist using homogenous proximity assay method. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

- Knaus *et al* (2007) Alpha2-adrenoceptor subtypes--unexpected functions for receptors and ligands derived from gene-targeted mouse models. *Neurochem Int* 51:277-281.
- Philipp *et al* (2002) Physiological significance of alpha(2)-adrenergic receptor subtype diversity: one receptor is not enough. *Am J Physiol Regul Integr Comp Physiol* 283:R287-R295.

FOR RESEARCH USE ONLY.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.