

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT α2C RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1436-1

Lot Number: C1436-1-101209

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

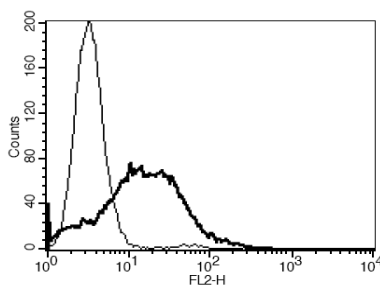
Transfection: Expression vector containing full-length human ADRA2C cDNA (GenBank Accession Number AY455666) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 µg/mL puromycin

Stability: Stable in culture for minimum of two months

Figure 3



Background: The human adrenergic α2c receptor is a 458-amino acid, 7-transmembrane protein. It belongs to the group of nine adrenoceptors that mediate the biological actions of the endogenous catecholamines adrenaline and noradrenaline. All three alpha2-receptor subtypes may operate as presynaptic inhibitory feedback receptors to control the release of neurotransmitters. The α2c receptor is of particular importance as it appears to regulate venous vasoconstriction.

Application: Functional assays

Figure 1

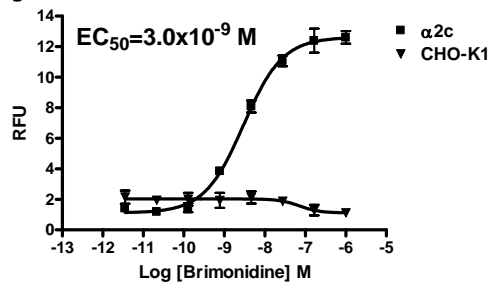


Figure 2

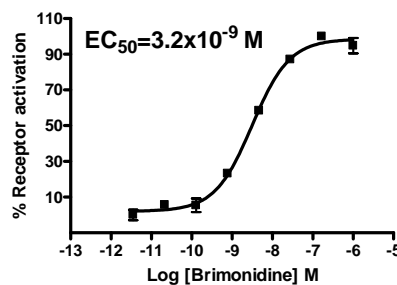


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Gyires *et al.* (2009) alpha(2)-Adrenoceptor subtypes-mediated physiological, pharmacological actions. *Neurochem Int* 55:447-453.

Knaus *et al* (2007) Alpha2-adrenoceptor subtypes--unexpected functions for receptors and ligands derived from gene-targeted mouse models. *Neurochem Int* 51:277-281.

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.