

MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT α2C RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1436-1

Lot Number: DC1436-1-072419

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cell Banker 2

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

Background: The human adrenergic α_{2c} 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 α_{2c} receptor is of particular importance as it appears to regulate venous vasoconstriction.

Application: Functional assays

Figure 1

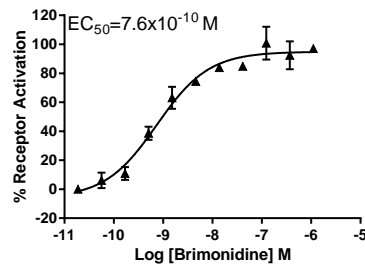


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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.

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