

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT M2 RECEPTOR

PRODUCT INFORMATION

Catalog Number: DC1023-1

Lot Number: DC1023-1-021721

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2 (Amsbio 11891)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human CHRM2 cDNA (GenBank Accession Number NM_001006632.1) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Data sheet

Background: The muscarinic M2 receptor is a 466-amino acid, 7-transmembrane protein. In heart muscle, M2 receptors represent the prevailing subtype of muscarinic receptors. They can also be found in neurons of central and peripheral nervous system. In neuronal cells, mainly on synaptic terminals, stimulation of the M2 autoreceptors is responsible for presynaptic muscarinic autoinhibition of acetylcholine release in both central and peripheral cholinergic neurons. Loss of function of these M2 receptors, as occurs in some patients with asthma and in animal models of inflammation, leads to an increase in vagally mediated hyperreactivity.

Application: Functional assays

Figure 1

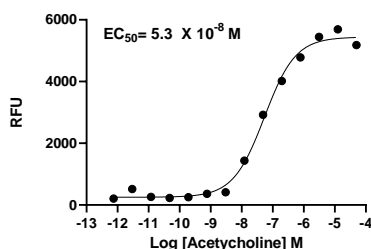


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Eglen (2006) Muscarinic receptor subtypes in neuronal and non-neuronal cholinergic function. *Auton Atacoid Pharmacol* 26:219-233.

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.