

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT M4 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1025-1

Lot Number: C1025-1-101209

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

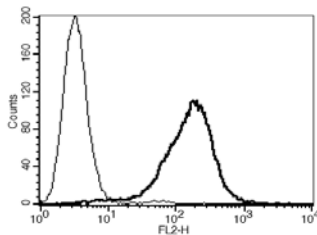
Transfection: Expression vector containing full-length human CHRM4 cDNA (GenBank accession number NM_000741.2) with FLAG tag sequence at N-terminus

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 muscarinic M4 receptor is a 479-amino acid, 7-transmembrane protein. In heart muscle, M4 receptors are coupled to specific K⁺ currents that regulate cardiac muscle contractions. They can also be found in tissues of the central nervous system and bladder.

Application: Functional assays
Figure 1

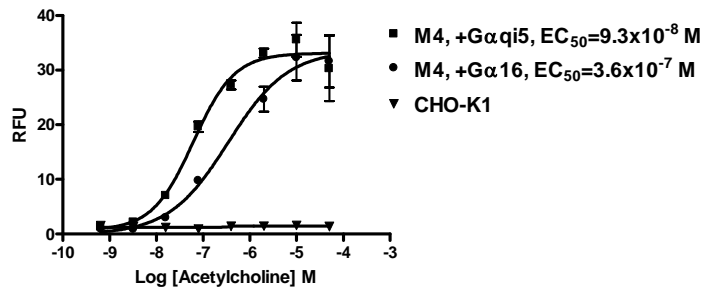


Figure 2

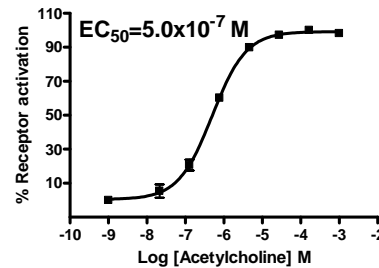


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent stimulation of intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 3.** Receptor expression on cell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

References:

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

Yeh *et al.* (2007) Atrial tachycardia induces remodeling of muscarinic receptors and their coupled potassium currents in canine left atrial and pulmonary vein cardiomyocytes. *Br J Pharmacol* 152:981-982.

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.