

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT M4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1025-1

Lot Number: C1025-1-07122013

Quantity: 10.65mg/ml, 1mg

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

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.

Figure 1

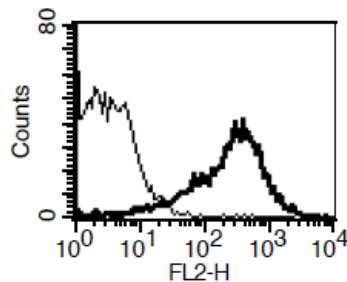


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Eglen (2006) Muscarinic receptor subtypes in neuronal and non-neuronal cholinergic function. *Auton Toxicol 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.