

**MULTISCREEN™ MEMBRANE PREPARATION
HUMAN RECOMBINANT μ OPIOID RECEPTOR**

PRODUCT INFORMATION

Catalog Number: MC1350-1a

Lot Number: MC1350-1a-07262012

Quantity: 1 vial (2.5 mg/ml)

Host cell: CHO-K1

Transfection: Expression vector containing full-length human OPRM1 cDNA (GenBank Accession Number: NM_000914.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen

Data sheet

Background: μ opioid receptor (MOR) is a G protein-coupled receptor for β-endorphin. The receptor activation inhibits neurotransmitter release by reducing calcium currents and increasing potassium conductance. MOR mediates positive reinforcement following direct (morphine) or indirect (alcohol, cannabinoids, nicotine) activation. MOR plays a genetic role in the control of gut inflammation. MOR-deficient mice are highly susceptible to colon inflammation, with a 50% mortality rate occurring 3 days after administration of TNBS that induces inflammation. MOR agonists regulate cytokine production and T cell proliferation and might be new therapeutic molecules in inflammatory bowel disease.

Application: [³⁵S] GTPγS Binding Assay

Figure 1

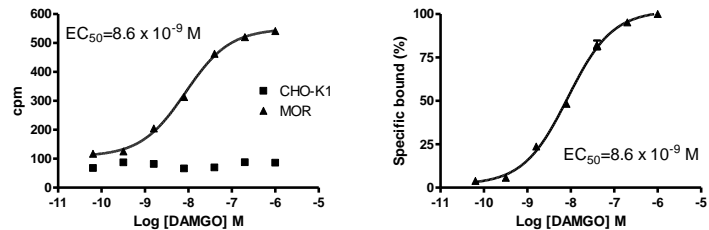


Figure 2

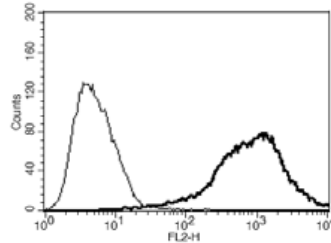


Figure 1. Dose-dependent [³⁵S]GTPγS Binding to OPRM1 membrane upon treatment with agonist using homogenous proximity assay method. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Chen *et al.* (1993) Molecular cloning and functional expression of a mu-opioid receptor from rat brain. *Mol Pharmacol* 44:8-12.

Contet *et al.* (2004) Mu opioid receptor: a gateway to drug addiction. *Curr Opin Neurobiol* 14:370-378.

Philippe *et al.* (2003) Anti-inflammatory properties of the mu opioid receptor support its use in the treatment of colon inflammation. *J Clin Invest* 111:1329-1338.

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.