

MULTISCREEN™ STABLE CELL LINE RAT RECOMBINANT MCHR1 RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1031-4

Lot Number: C1031-4-092410

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gαq5

Transfection: Full-length rat Mchr1 cDNA (GenBank Accession Number NM_031758)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 100 µg/mL zeocin, 100 µg/mL G418

Stability: Stable in culture for minimum of two months

Data sheet

Background: MCHR1 or GPR24 is a receptor for melanin-concentrating hormone (MCH). MCH-1R is one of the two MCH receptors identified in human. The primary function of both MCH receptors, MCHR1 and MCHR2, is to regulate skin color. In addition, they also regulate food intake, energy expenditure, behavior and emotion. Blockade of MCHR1 could provide a viable approach for treatment of obesity and certain neurological disorders.

Application: Functional assays and binding assays

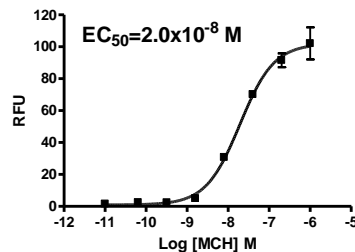


Figure legend: Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Mori *et al.* (2001) Cloning of a novel G protein-coupled receptor, SLT, a subtype of the melanin-concentrating hormone receptor. *Biochem Biophys Res Commun* 283:1013-1018.

Shi (2004) Beyond skin color: emerging roles of melanin-concentrating hormone in energy homeostasis and other physiological functions. *Peptides* 25:1605-1611.

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. June 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.