

**MULTISCREEN™**  
**HUMAN ENDOGENOUS MCH1 (GPR24) RECEPTOR**

---

**PRODUCT INFORMATION**

**Catalog Number:** C1031-1

**Lot Number:** C1031-1-061505

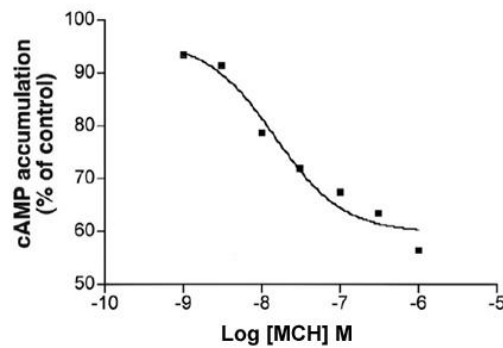
**Cell:** YS35 naturally expressing human MCH1R

**Propagation Medium:** RPMI1640, 10% FBS

**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



**Figure legend:** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand.

**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 SCREENING BY MULTISPAN INC. ONLY.**  
**NOT SOLD AS PRODUCT.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

26219 Eden Landing Road, Hayward, California 94545-3718 Phone: (510) 887-0817 [info@multispaninc.com](mailto:info@multispaninc.com)