

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT MCH2 (GPR145) RECEPTOR**

---

**PRODUCT INFORMATION**

**Catalog Number:** C1032

**Lot Number:** C1032-061505

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO dhfr<sup>-</sup>

**Transfection:** Full-length human MCH-2R/GPR145 cDNA (GenBank Accession Number NM\_032503.1)

**Recommended Storage:** Liquid nitrogen upon receiving

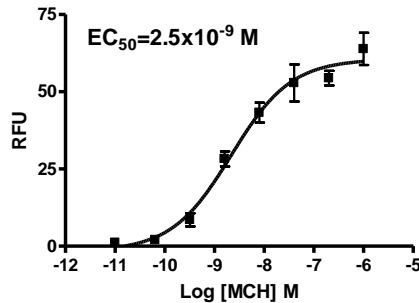
**Propagation Medium:** Alpha-MEM, 10% FBS, 400 µg/mL G418

**Stability:** Stable in culture for minimum of two months

**Data sheet**

**Background:** Melanin-concentrating hormone receptor 2 (MCHR2, or GPR145) is the second MCH receptor 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.

**Application:** Functional 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.