

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT MCH1 RECEPTOR (GPR24)**

---

**PRODUCT INFORMATION**

**Catalog Number:** C1031-3

**Lot Number:** C1031-3-073005

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

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

**Host cell:** HEK293T Gaqi5

**Transfection:** Full-length human MCHR1 cDNA (GenBank Accession Number NM\_005297)

**Recommended Storage:** Liquid nitrogen upon receiving

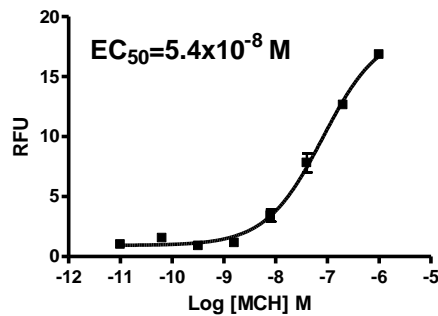
**Propagation Medium:** DMEM, 10% FBS, 200 µg/mL hygromycin, 100 µg/mL zeocin

**Stability:** Stable after minimum of two months continuous growth

**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 stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

**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.**

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