

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

PRODUCT INFORMATION

Catalog Number: C1032

Lot Number: C1032-073005

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO dhfr⁻

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

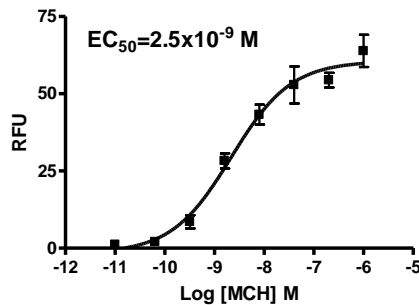


Figure legend: Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.