

PRODUCT DATASHEET
ChemiScreen™ MC₁ Melanocortin Membrane Preparation

CATALOG NUMBER:	HTS156M	QUANTITY:	200 units
LOT NUMBER:	JH1909399	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL

BACKGROUND:

The melanocortin-1 receptor (MC₁) is a member of the G-protein-coupled receptor family of proteins that is responsible for regulating skin and hair color. The MC₁ receptor is regulated by the physiological agonists alpha-melanocyte-stimulating hormone (alphaMSH) and adrenocorticotrophic hormone (ACTH), and antagonist agouti signaling protein (ASP) (Rouzud *et al.*, 2005). In 1995 a landmark study demonstrated that over 80% of humans with red hair or fair skin have a dysfunctional variant of the MC₁ gene (Valverde *et al.*, 1995). MC₁ is, therefore, a major determinant of sun sensitivity and genetic risk factor for melanoma and non-melanoma skin cancer. MC₁ membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level of GPCR surface expression; thus, they are ideal HTS tools for screening of agonists and antagonists of MC₁. The membrane preparations exhibit a K_d of 1.3nM for [¹²⁵I]-[Nle⁴, D-Phe⁷]-α-MSH (NDP-αMSH). With 0.3 nM [¹²⁵I]-NDP-αMSH, 5μg/well MC₁ Membrane Prep typically yields greater than 50-fold signal-to-background ratio.

APPLICATIONS:

Radioligand binding assay and GTPγS binding

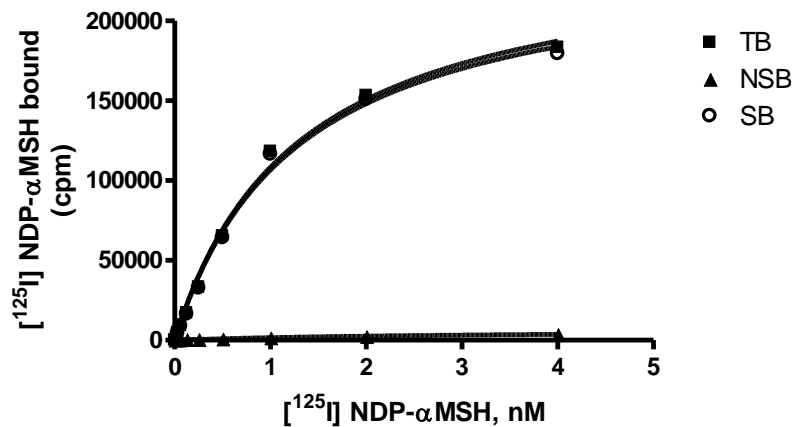


Figure 1. Saturation binding for MC₁. 5 μg/well MC₁ Membrane Preparation was incubated with increasing amount of [¹²⁵I]-NDP-αMSH in the absence (total binding, TB) or presence (nonspecific binding, NSB) of greater than 500-fold excess unlabeled NDP-αMSH. Specific binding (SB) was determined by subtracting NSB from TB.

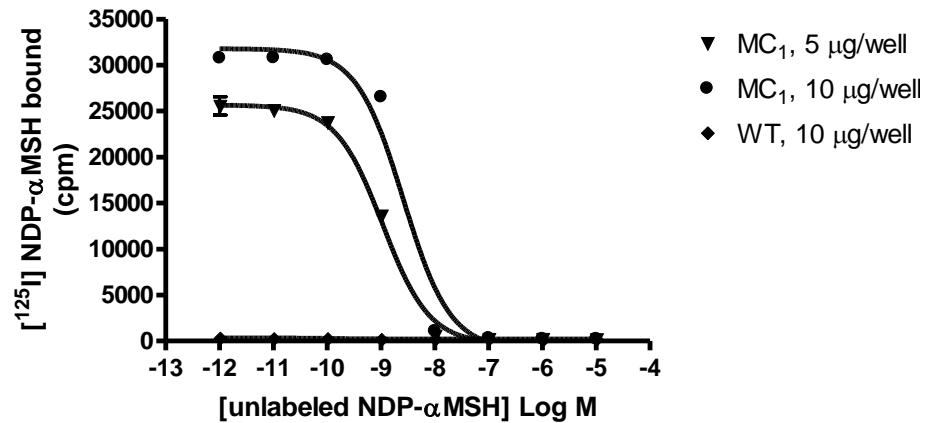


Figure 2. Competition binding for MC₁. MC₁ Membrane Preparation (5 or 10 μg/well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.3 nM [¹²⁵I]-NDP-αMSH and increasing concentrations of unlabeled NDP-αMSH, and more than 50- fold signal:background was obtained.

SPECIFICATIONS: 1 unit = 5 μg
 B_{max} for [¹²⁵I]-NDP-αMSH binding: 13.7 pmol/mg protein
 K_d for [¹²⁵I]-NDP-αMSH binding: ~1.3 nM
 Signal:background: >50-fold

TRANSFECTION: Full-length human MC1R cDNA encoding MC₁ (Accession Number: NM_002386).

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous MC₁ expression.

RECOMMENDED ASSAY CONDITIONS: Membranes are mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a nonbinding 96-well plate, and incubated for 1-2 h. Prior to filtration, an FC 96-well harvest plate (EMD Millipore cat. # MAHF C1H) is coated with 0.33% polyethyleneimine for 30 min, then washed with 50mM HEPES, pH 7.4, 500mM NaCl. Binding reaction is transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The plate is dried and counted.

Binding buffer: 50 mM Hepes, pH 7.4, 5 mM MgCl₂, 1 mM CaCl₂, filtered and stored at 4°C

Radioligand: [¹²⁵I]-NDP-αMSH (Perkin Elmer#: NEX-352)

Wash Buffer: 50 mM Hepes, pH 7.4, 500mM NaCl, filtered and stored at 4°C.

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 50-fold signal:background with ¹²⁵I labeled NDP-αMSH at 0.3 nM.

PRESENTATION: Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.
 Packaging method: Membranes protein were adjusted to 1 mg/mL in packaging buffer, and dispensed at 1 mL/vial. Vials were rapidly frozen, and stored at -80°C.

STORAGE/HANDLING:

Store at -70°C . Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

REFERENCES:

1. Valverde P *et al.* (1995) Variants of the melanocyte-stimulating hormone receptor gene are associated with red hair and fair skin in humans. *Nat. Genet.* 11: 328-30.
2. Rouzaud F *et al.* (2005) MC1R and the response of melanocytes to ultraviolet radiation. *Mutat. Res.* 571: 133-152.

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services