

PRODUCT DATASHEET

ChemiScreen™ MC₅ Melanocortin Membrane Preparation

CATALOG NUMBER:	HTS155M	QUANTITY:	200 units
LOT NUMBER:	SC20190807	VOLUME/CONCENTRATION:	1 mL, 2 mg/mL

BACKGROUND: The melanocortin system consists of five seven-transmembrane spanning G-protein coupled receptors (MC₁-MC₅). Among the five members of the melanocortin receptor family, MC₂ and MC₅ are expressed in peripheral tissues. The MC₂ receptor (ACTH receptor) is almost exclusively expressed in the adrenal cortex whereas MC₅ is expressed in several organs including the adrenal cortex. Studies have shown that targeted disruption of the MC5R gene produced mice with a severe defect in water repulsion and thermoregulation caused by decreased production of sebaceous lipids (Chen *et al.*, 1997). Data show a requirement for the MC₅ in multiple exocrine glands for the production of numerous products, indicative of a coordinated system for regulation of exocrine gland function by melanocortin peptides. (Entwistle *et al.*1990). MC₅ 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 MC₅.

APPLICATIONS: Radioligand Binding Assay

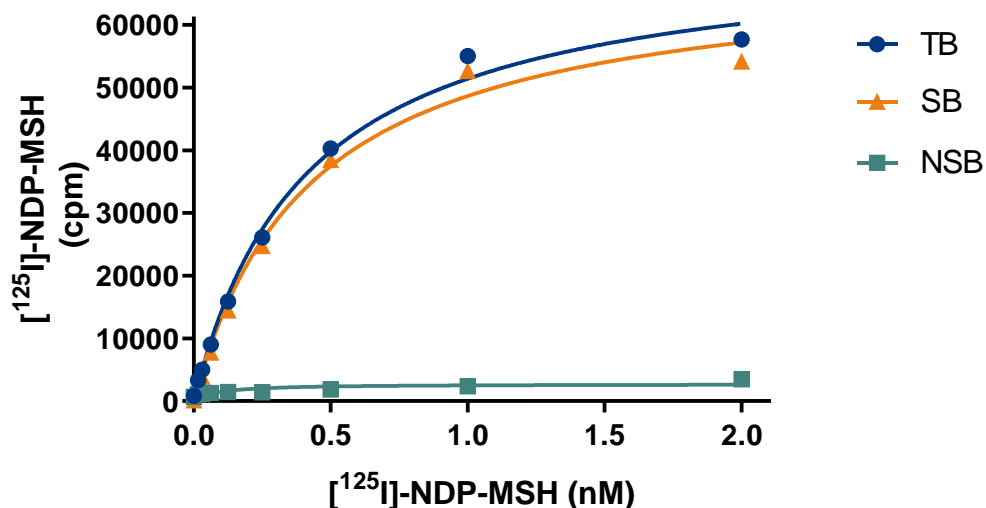


Figure 1. Saturation Binding for MC₅. 10 μg/well MC₅ Membrane Preparations were incubated with increasing amount of [¹²⁵I]-NDP-αMSH in the absence (total binding, TB) or presence (nonspecific binding, NSB) of greater than 500-fold excess unlabeled NDP-αMSH. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

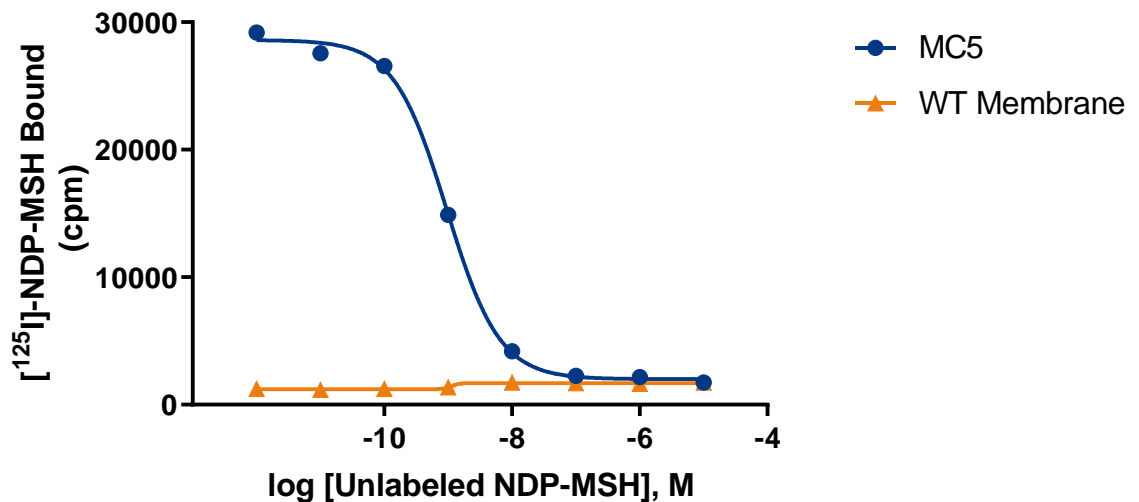


Figure 2. Competition Binding for MC₅. MC₅ Membrane Preparation and Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1), each at 10 µg/well, were incubated with 0.3 nM [¹²⁵I]-NDP-αMSH and increasing concentrations of unlabeled NDP-αMSH, and more than 15- fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max} for [¹²⁵I]-NDP-αMSH binding: 2.84 pmol/mg
 K_d for [¹²⁵I]-NDP-αMSH binding: 0.43 nM
 Signal:background: ≥15-fold

TRANSFECTION: Full-length human MC5R cDNA encoding MC₅ (Accession Number: NM_005913).

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous MC₅ 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 50 mM HEPES, pH 7.4, 500 mM 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 (PerkinElmer#: NEX-352)

Wash Buffer: 50 mM HEPES, pH 7.4, 500 mM 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] 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: Membrane proteins were adjusted to the indicated concentration in 1 ml packaging buffer, 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. Avoid repeated freeze/thaw cycles.

REFERENCES:

1. Chen W *et al.* (1997). Exocrine gland dysfunction in MC5-R-deficient mice: evidence for coordinated regulation of exocrine gland function by melanocortin peptides. *Cell* 91:789-798.
2. Entwistle ML *et al.* (1990). Characterization of functional melanotropin receptors in lacrimal glands of the rat. *Peptides* 11:477-483.

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.