

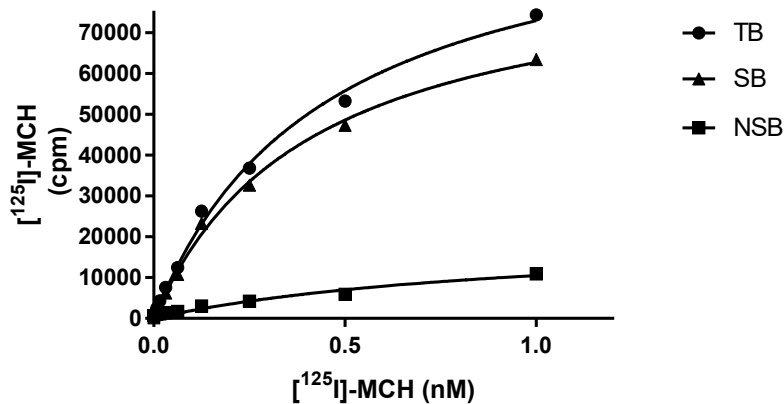
PRODUCT DATASHEET

ChemiScreen™ MCH1 Membrane Preparation

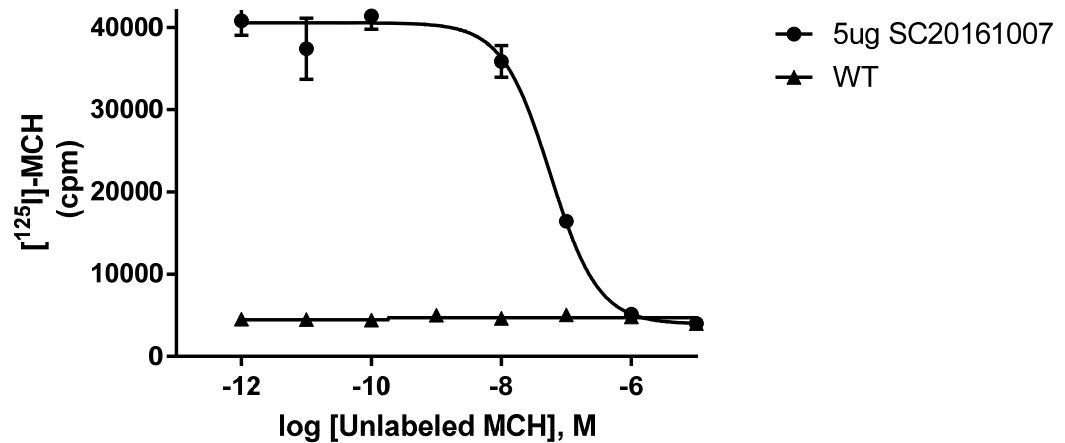
**CATALOG NUMBER:** HTS025M                      **QUANTITY:** 200 units  
**LOT NUMBER:** SC20161007                      **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

**BACKGROUND:** Melanin-concentrating hormone (MCH) is a cyclic, 19 amino acid peptide produced in the lateral hypothalamus and other regions of the mammalian brain (Bittencourt *et al.*, 1992). When administered intracerebrally or overexpressed transgenically in rodents, MCH stimulates food intake and weight gain (Ludwig *et al.*, 2001). A receptor for MCH, previously identified as an orphan receptor SLC-1 and renamed MCH<sub>1</sub>, has been identified in humans and rodents. Consistent with a role in hypothalamic control of feeding behavior, MCH<sub>1</sub> is expressed in in the ventromedial and dorsomedial nuclei of the hypothalamus (Chambers *et al.*, 1999). Genetic ablation and pharmacological antagonism of MCH<sub>1</sub> in rodents results in resistance to obesity caused by a high fat diet or leptin deficiency (Borowsky *et al.*, 2002; Segal-Lieberman, *et al.*, 2003; Shearman *et al.*, 2003; Marsh *et al.*, 2002; Chen *et al.*, 2002). Thus, MCH<sub>1</sub> is an attractive target for obesity. MCH<sub>1</sub> membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level GPCR surface expression; thus, they are ideal HTS tools for screening for agonists and antagonists of MCH<sub>1</sub>.

**APPLICATIONS:** Radioligand binding assay



**Figure 1. Saturation binding for MCH<sub>1</sub>.** 5.0 ug/well MCH<sub>1</sub> Membrane Preparation was incubated with increasing amount of [<sup>125</sup>I]-MCH in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled MCH. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for MCH<sub>1</sub>.** 5 ug/well MCH<sub>1</sub> Membrane Preparation and Wild-Type Chem-1 membrane preparation (10 μg/well; catalog # HTS000MC1) were incubated with 0.4 nM [<sup>125</sup>I]-MCH and increasing concentrations of unlabeled MCH, and more than 8-fold signal:background was obtained. Representative sample data.

**SPECIFICATIONS:** 1 unit = 5 μg  
 B<sub>max</sub>: 195.96 pmol/mg protein  
 K<sub>d</sub>: ~0.41 nM  
 Signal:background: >10-fold

**TRANSFECTION:** Full length human MCHR1 cDNA encoding MCH<sub>1</sub> (Accession number NM\_005297)

**Species:** Human

**HOST CELLS:** Chem-1, an adherent mammalian cell line with no detectable endogenous MCH1 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, 0.5% BSA. 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 MgCl<sub>2</sub>, 1 mM CaCl<sub>2</sub>, 0.2% BSA, filtered and stored at 4°C

**Radioligand:** [<sup>125</sup>I]-MCH (Perkin Elmer CUSTOM)

**Wash Buffer:** 50 mM Hepes, pH 7.4, 500mM NaCl, 0.1% BSA, filtered and stored at 4°C

One package contains enough membranes for at least 200 assays (units), where an unit is the amount of membrane that will yield greater than 8-fold signal:background with <sup>125</sup>I-labeled MCH at 0.4 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. Do not freeze and thaw.
- REFERENCES:**
1. Bittencourt J.C., *et al.* (1992) The melanin-concentrating hormone system of the rat brain: an immuno- and hybridization histochemical characterization. *J. Comp. Neurol.* 319: 218-45.
  2. Borowsky B., *et al.* (2002) Antidepressant, anxiolytic and anorectic effects of a melanin-concentrating hormone-1 receptor antagonist. *Nat. Med.* 8: 825-30.
  3. Chambers J., *et al.* (1999) Melanin-concentrating hormone is the cognate ligand for the orphan G-protein-coupled receptor SLC-1. *Nature* 400: 261-5.
  4. Chen Y., *et al.* (2002) Targeted disruption of the melanin-concentrating hormone receptor-1 results in hyperphagia and resistance to diet-induced obesity. *Endocrinology* 143: 2469-77.
  5. Ludwig D.S., *et al.* (2001) Melanin-concentrating hormone overexpression in transgenic mice leads to obesity and insulin resistance. *J. Clin. Invest.* 107: 379-86
  6. Marsh D.J., *et al.* (2002) Melanin-concentrating hormone 1 receptor-deficient mice are lean, hyperactive, and hyperphagic and have altered metabolism. *Proc. Natl. Acad. Sci. USA* 99: 3240-5.
  7. Segal-Lieberman G., *et al.* (2003) Melanin-concentrating hormone is a critical mediator of the leptin-deficient phenotype. *Proc. Natl. Acad. Sci. USA* 100: 10085-90.
  8. Shearman L.P., *et al.* (2003) Chronic MCH-1 receptor modulation alters appetite, body weight and adiposity in rats. *Eur. J. Pharmacol.* 475: 37-47.

**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