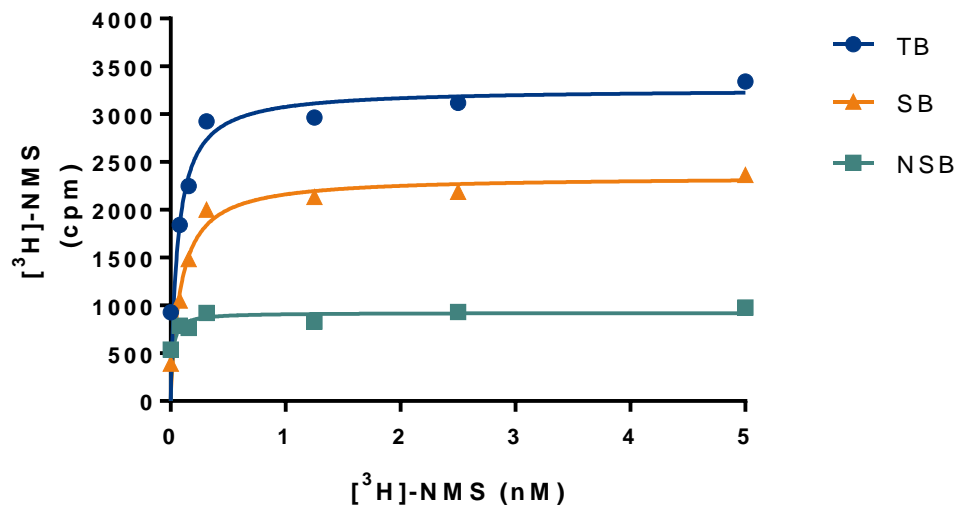


**PRODUCT DATASHEET**
**ChemiScreen™ M<sub>4</sub> Muscarinic Acetylcholine Membrane Preparation**

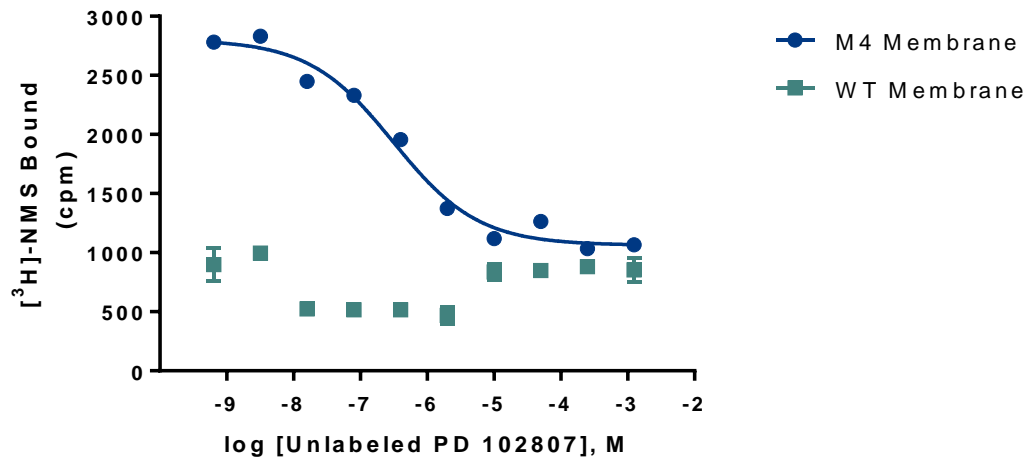
**CATALOG NUMBER:** HTS117M                      **QUANTITY:** 200 units  
**LOT NUMBER:** SC20181211                      **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

**BACKGROUND:** The muscarinic acetylcholine receptor (mAChR) family consists of five GPCRs that mediate some of the neurotransmission functions of acetylcholine in the CNS and the periphery. The M<sub>1</sub>, M<sub>3</sub> and M<sub>5</sub> receptors couple to Gq to mobilize intracellular calcium, whereas the M<sub>2</sub> and M<sub>4</sub> receptors couple to Gi/o to inhibit cAMP production (Caulfield and Birdsall, 1998). Muscarinic M<sub>4</sub> receptor is known to be abundantly expressed in the striatum (Levey 1993). Consistent with this fact, locomotor activity was significantly increased in M<sub>4</sub> receptor-deficient mice (Gomez et al. 1999). M<sub>4</sub> 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 M<sub>4</sub>.

**APPLICATIONS:** Radioligand binding assay



**Figure 1. Saturation binding for M<sub>4</sub>.** 10 μg/well M<sub>4</sub> Membrane Preparation was incubated with increasing amount of <sup>3</sup>H-labeled NMS in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 1000-fold excess unlabeled PD 102807. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.



**Figure 2. Competition binding for M<sub>4</sub>.** 10 µg/well M<sub>4</sub> Membrane Preparation and 10 µg/well wild-type Chem-1 Membrane Preparation (catalog # HTS000MC1) were incubated in a 96-well plate with 0.4 nM <sup>3</sup>H-labeled NMS and increasing concentrations of unlabeled PD 102807. More than 2.5-fold signal:background was obtained. Representative sample data.

**SPECIFICATIONS:** 1 unit = 10 µg  
 $B_{max}$  for [<sup>3</sup>H]-NMS binding: 2.64 pmol/mg protein  
 $K_d$  for [<sup>3</sup>H]- NMS binding: ~0.09 nM  
Signal:background: >2.5-fold

**TRANSFECTION:** Human M4 cDNA (Accession Number: NM\_00741)

**Species:** Human

**HOST CELLS:** Chem-4, an adherent cell line expressing the promiscuous G-protein, Gα15.

**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, 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>3</sup>H]-NMS. (Perkin Elmer#:NET636 )

**Wash Buffer:** 50 mM HEPES, pH 7.4, 500 mM NaCl, 0.1% BSA, 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 2.5-fold signal:background with [<sup>3</sup>H]-NMS.

**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  $-70^{\circ}\text{C}$ . Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

**REFERENCES:**

1. Caulfield MP and Birdsall NJM (1998) International Union of Pharmacology. XVII. Classification of muscarinic acetylcholine receptors. *Pharmacol. Rev.* 50: 279-290.
2. Levey AI (1993). Immunological localization of M1-M5 muscarinic acetylcholine receptors in peripheral tissue and brain. *Life Sci.* 52: 441-448.
3. Gomeza J et al. (1999) Enhancement of D1 dopamine receptor-mediated locomotor stimulation in M4 muscarinic acetylcholine receptor knockout mice. *Proc. Natl. Acad. Sci. USA* 96: 10483-10488.

**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