

PRODUCT DATASHEET
ChemiScreen™ M₂ Muscarinic Acetylcholine Membrane Preparation

CATALOG NUMBER:	HTS115M	QUANTITY:	200 units
LOT NUMBER:	SC20180621	VOLUME/CONCENTRATION:	1 mL, 2.0 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 M₁, M₃ and M₅ receptors couple to G_q to mobilize intracellular calcium, whereas the M₂ and M₄ receptors couple to G_{i/o} to inhibit cAMP production (Caulfield and Birdsall, 1998). In urinary bladder trachea and stomach, M₂ augments the function of M₃ in promoting contractility, and activation of M₂ serves to counteract relaxation induced by increased cAMP levels (Ehlert et al., 2005; Wess, 2004). In addition, the ability of mAChR agonists to decrease heart rate appears to be mediated primarily by M₂. Agonists of mAChRs induce tremor, hypothermia, corticosterone release, and analgesia; each of these functions is mediated at least in part by M₂ (Wess, 2004). The M₂ 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 M₂ interactions with its ligands. The membrane preparations exhibit a K_d of 0.76 nM for [³H]-Scopolamine methyl chloride (NMS). With 10 μg/well M₂ Membrane Prep and 0.5 nM [³H]-NMS, greater than 3-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand binding assay

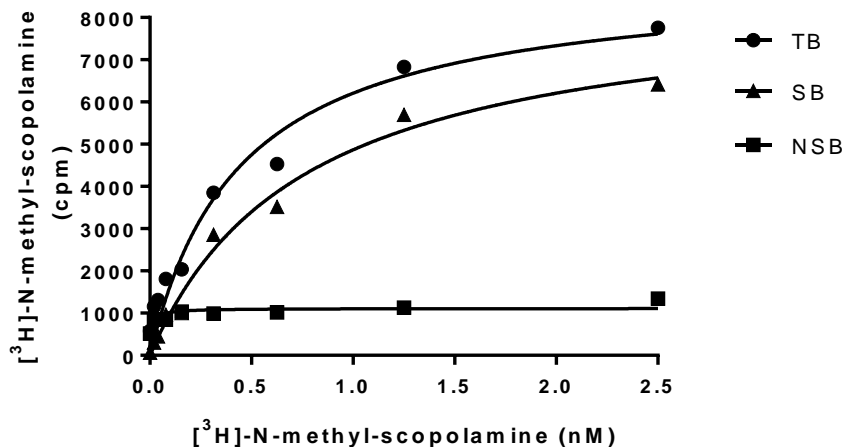


Figure 1. Saturation binding for M₂. 10 μg/well M₂ Membrane Preparation was incubated with increasing amount of ³H-labeled NMS in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled Atropine. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

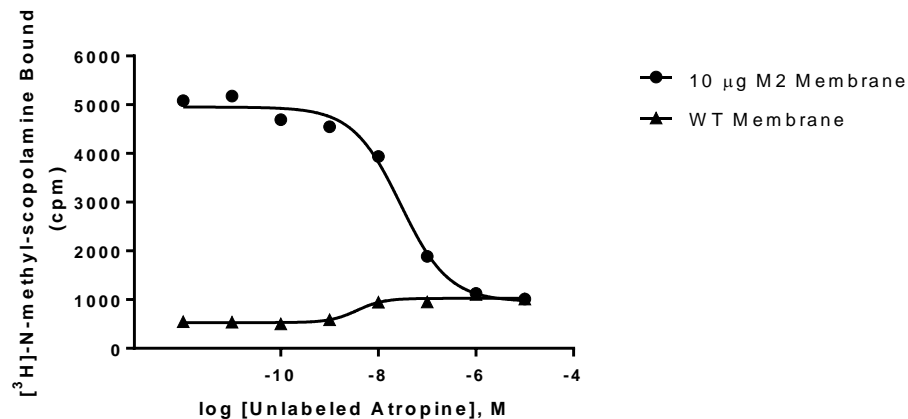


Figure 2. Competition binding for M₂: 10 µg/well M₂ Membrane Preparation (HTS115M) or 10 µg/well Chem-1 wild-type membrane (HTS000MC1) were incubated with 0.5 nM ³H-labeled NMS and increasing concentrations of unlabeled Atropine. More than 3- fold signal:background ratio was obtained. Sample data from a representative lot.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max} for [³H]- N-methyl-scopolamine binding: 9.65 pmol/mg protein
 K_d for [³H]- N-methyl-scopolamine binding: ~0.76 nM
 Signal:Background: >3-fold

Species: Full-length human CHRM2 cDNA encoding M₂ (Accession Number: NM_000739)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous M₂ 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 4mM Na₂HPO₄, 1mM KH₂PO₄, pH7.4. 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: 4mM Na₂HPO₄, 1mM KH₂PO₄, pH7.4, filtered and stored at 4°C.

Radioligand: [³H] NMS (Perkin Elmer#:NEX-636)

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 3-fold signal:background with ³H-labeled N-methyl-scopolamine at 0.5 nM.

PRESENTATION:

Liquid in packaging buffer: 50 mM Tris, pH 7.4, 10% glycerol and 1% BSA with no preservatives.

Packaging method: Membranes protein was adjusted to the indicated concentration in 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. Caulfield MP and Birdsall NJM (1998) International Union of Pharmacology. XVII. Classification of muscarinic acetylcholine receptors. *Pharmacol. Rev.* 50: 279-290.
2. Ehlert FJ *et al.* (2005) The M₂ muscarinic receptor mediates contraction through indirect mechanisms in mouse urinary bladder. *J. Pharmacol. Exp. Ther.* 313: 368-378.
3. Wess J (2004) Muscarinic acetylcholine knockout mice: novel phenotypes and clinical implications. *Annu. Rev. Pharmacol. Toxicol.* 44: 423-450.

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