

PRODUCT DATASHEET
ChemiScreen™ M₃ Muscarinic Acetylcholine Membrane Preparation

CATALOG NUMBER:	HTS116M	QUANTITY:	200 units
LOT NUMBER:	SC20180830	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 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). M₃ is expressed prominently in smooth muscle, and plays a primary role in mediating mAChR agonist-induced contractility. Mice lacking M₃ have dilated pupils, which indicates a role for M₃ in regulating tone of the pupillary sphincter muscle. In addition, M₃ plays a role in feeding, as indicated by the lean and hypophagic phenotype of M₃-null mice (Wess, 2004). M₃ 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 antagonists of M₃ interactions with 4-DAMP. The membrane preparations exhibit a K_d of 0.49 nM for [³H]-4-DAMP. With 10 μg/well M₃ Membrane Prep and 0.75 nM [³H]-4-DAMP, a greater than 2.5-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand binding assay

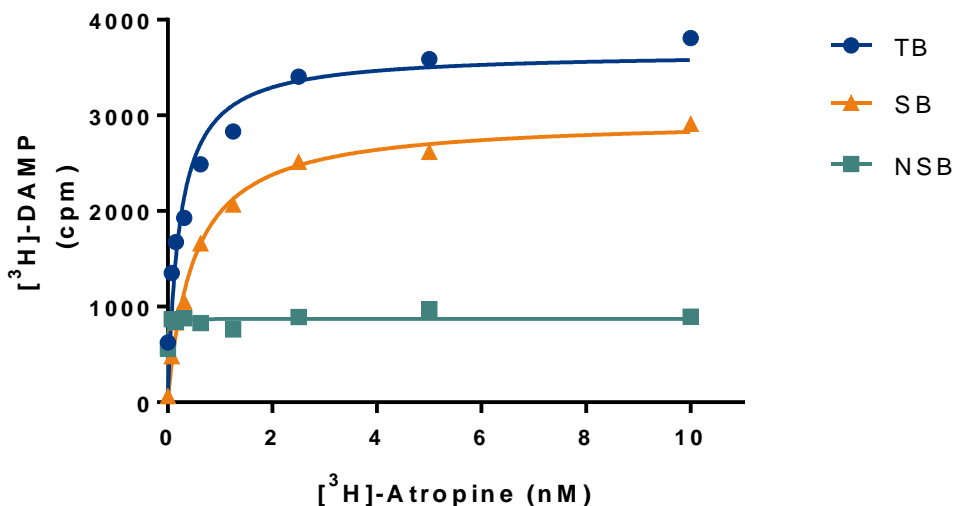


Figure 1. Saturation binding for M₃. 10 μg/well M₃ Membrane Preparation was incubated with increasing amount of [³H]-4-DAMP in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 1000-fold excess unlabeled atropine. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

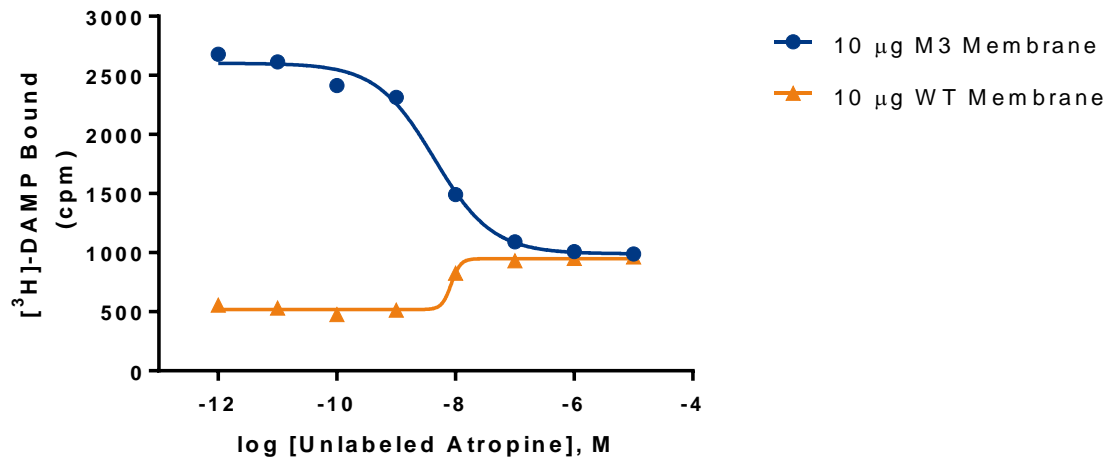


Figure 2. Competition binding for M₃. M₃ Membrane Preparation (10 µg/well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.75 nM [³H]-4-DAMP and increasing concentrations of unlabeled atropine, and more than 2.5- fold signal:background was obtained. Sample data from a representative lot.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max}: 3.22 pmol/mg protein
 K_d: ~0.49 nM
 Signal:background: >2.5-fold

TRANSFECTION: Full-length human CHRM3 cDNA encoding M₃ (Accession Number: NM_000740)

Species: Human

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous M₃ 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, a GF/C 96-well filter plate 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 MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C

Radioligand: [³H] 4-DAMP (Perkin Elmer # NET1040)

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 a unit is the amount of membrane that will yield greater than 2.5-fold signal:background with ³H-labeled 4-DAMP.

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. Caulfield MP and Birdsall NJM (1998) International Union of Pharmacology. XVII. Classification of muscarinic acetylcholine receptors. *Pharmacol. Rev.* 50: 279-290.
 2. 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