

PRODUCT DATASHEET
ChemiScreen™ H₁ HISTAMINE Membrane Preparation

CATALOG NUMBER: HTS050M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

BACKGROUND: Histamine exerts its biological effects through a family of four GPCRs. Stimulation of the H₁ histamine receptor causes smooth muscle contraction, induces vascular permeability, and stimulates catecholamine release from the adrenal medulla. H₁ is also widely expressed in the brain, and activation of H₁ excites several types of neurons by blocking potassium conductance (Hill *et al.*, 1997). H₁ 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 H₁ interactions and its ligands. The membrane preparations exhibit a K_d of 4.5 nM for [³H]-Pyrilamine. With 3 nM [³H]-Pyrilamine, 10 µg/well of H₁ Membrane Prep typically yields a greater than 4-fold signal-to-background ratio.

APPLICATIONS: Radioligand Binding Assay

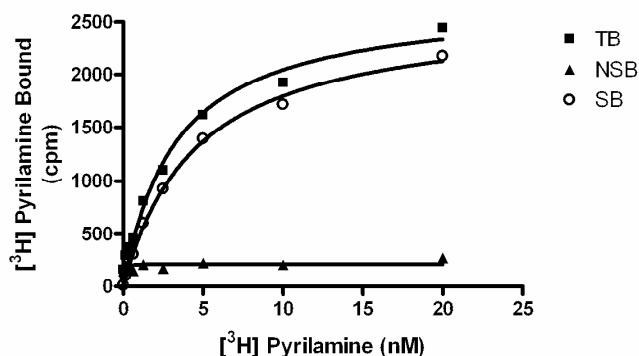


Figure 1. Saturation Binding for H₁. 10 µg/well of H₁ Membrane Preparation were incubated with increasing amounts of [³H]-Pyrilamine in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled Pyrilamine. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample of lot SC.

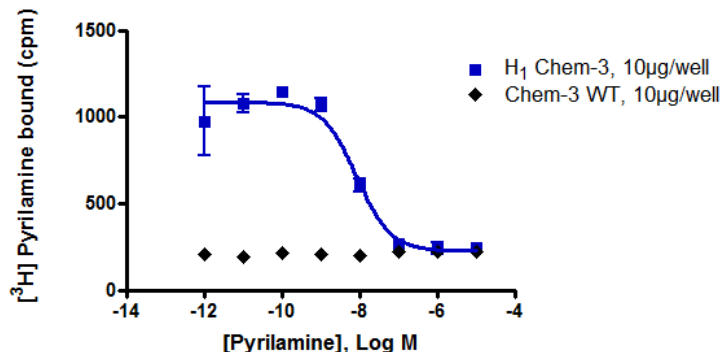


Figure 2. Competition Binding for H₁. 10 µg/well of H₁ or Wild Type Chem-3 (Catalog # HTS000MC3) Membrane Preparation were incubated with 3 nM [³H]-Pyrilamine and increasing concentrations of unlabeled Pyrilamine and were subjected to filtration binding. More than a 4-fold signal:background ratio was obtained. The data are from a representative sample of lot SC.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max} for [³H]-Pyrilamine Binding: 8.3 pmol/mg protein
 K_d for [³H]-Pyrilamine Binding: 4.5 nM
 Signal:Background: ≥4-fold

TRANSFECTION: Full-length human HRH1 cDNA encoding H₁ (Accession Number: NM_000861)

HOST CELLS: Chem-3, a suspension mammalian cell line without any endogenous H₁ expression.

RECOMMENDED ASSAY CONDITIONS: Membranes were mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a non-binding 96-well plate, and incubated for 2 h at room temperature. Prior to filtration, an FC 96-well harvest plate was coated with 0.33% polyethyleneimine for 30 min, then washed with 50 mM HEPES, pH 7.4, 0.5% BSA. The binding reactions were transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The wells were then dried and counted for determination of receptor-associated radioligand binding.

Binding Buffer: 50 mM Tris, pH 7.4, 10 mM MgCl₂, 1 mM EDTA, filtered and stored at 4°C.

Radioligand: [³H]-Pyrilamine (PerkinElmer#: NET594)

Wash Buffer: 50 mM Tris, pH 7.4, 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 a 4-fold signal:background ratio with [³H]-Pyrilamine.

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

Membranes proteins were adjusted to 2 mg/mL in packaging buffer, dispensed at 1 mL per vial, 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. Hill SJ *et al.* (1997). International Union of Pharmacology. XIII. Classification of histamine receptors. *Pharmacol. Rev.* 49:253-278.

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.