

PRODUCT DATASHEET
ChemiScreen™ A₁ Adenosine Membrane Preparation

CATALOG NUMBER: HTS047M **QUANTITY:** 200 units
LOT NUMBER: SC20181219 **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

BACKGROUND: Extracellular adenosine mediates a multitude of biological effects, including wakefulness, antiarrhythmia, bronchoconstriction and response to ischemia and oxidative stress. A family of four GPCR adenosine receptors, A₁, A_{2A}, A_{2B} and A₃, is responsible for these effects. The A₁ receptor, which couples to Gi/o, is most highly expressed in brain, and mediates endogenous antinociception and neuronal response to hypoxia (Fredholm et al., 2001). A₁ is also expressed in kidney, where it contributes to tubuloglomerular feedback. A₁ 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 at A₁. The membrane preparations exhibit a K_d of 4.6 nM for [³H]-Cyclopentyl-1, 3-dipropylxanthine, 8-[dipropyl-2, 3-3H (N)]. With 10 µg/well A₁ Membrane Prep and 20 nM [³H]-DPCPX, a greater than 6-fold signal-to-background ratio is obtained.

APPLICATIONS: Radioligand binding assay

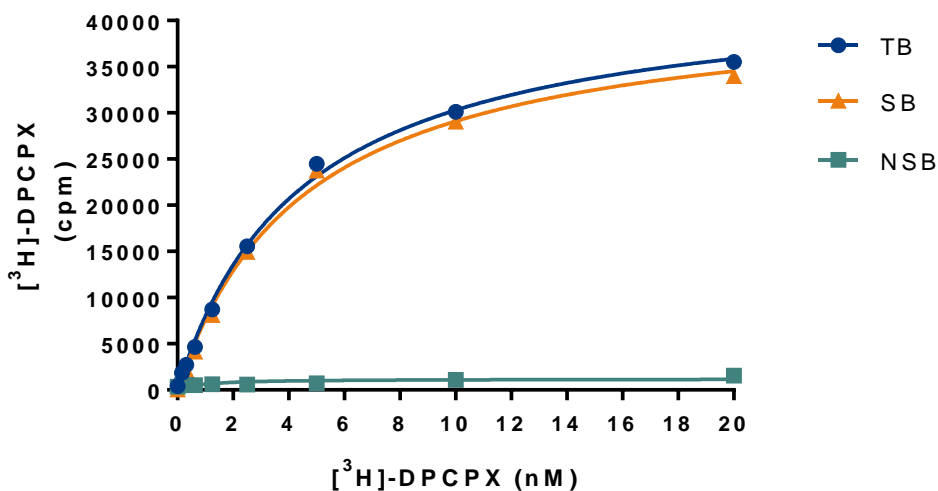


Figure 1. Saturation binding for A₁. 10 µg/well A₁ Membrane Preparation was incubated with increasing amount of ³H-labeled DPCPX in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled DPCPX. Specific binding (SB) was determined by subtracting NSB from TB.

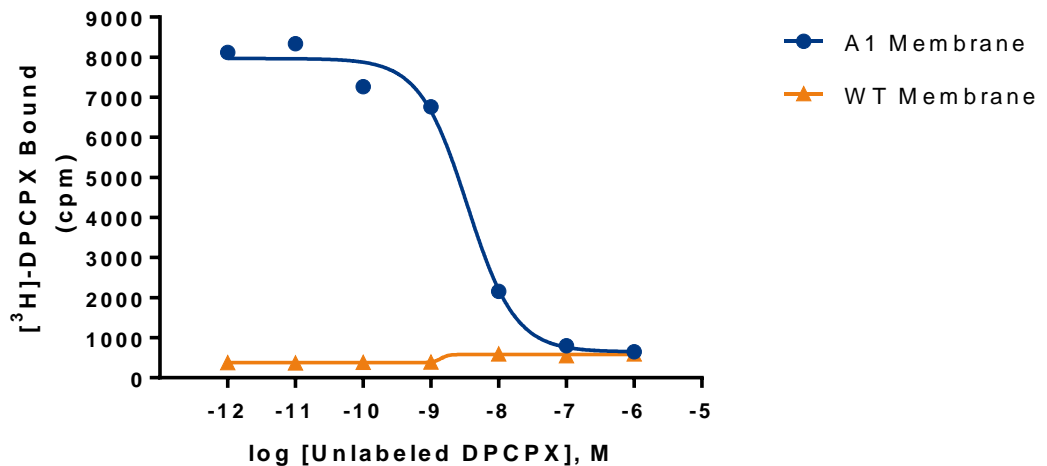


Figure 2. Competition binding for A₁. 10 µg/well A₁ Membrane Preparation (A₁; HTS047M) and wild type Membrane Preparation (WT) were incubated with 1 nM ³H-labeled DPCPX and increasing concentrations of unlabeled DPCPX. More than 6-fold signal:background was obtained with A₁ membranes, whereas no binding was observed with wild-type membranes.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max}: 23.28 pmol/mg
 K_d: 4.6 nM
 Signal:Background: >6-fold

Species: Human A₁ (Accession number S45235)

HOST CELLS: Chem-3, a suspension mammalian cell line without any endogenous A₁ 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 (Millipore cat. # MAHF C1H) is coated with 0.33% polyethyleneimine for 30 min, and then washed with 50mM Tris, pH 7.4. 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: 50 mM Tris, pH 7.4, 10 mM MgCl₂, 1 mM EDTA, filtered and stored at 4°C.

Radioligand: [³H] DPCPX (Perkin Elmer# :NET974)

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 6-fold signal:background with ³H-labeled DPCPX at 1 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 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. Fredholm, BB et al. (2001) International Union of Pharmacology. XXV. Nomenclature and classification of adenosine receptors. Pharmacol. Rev. 53: 527-552.

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