

PRODUCT DATASHEET
ChemiScreen™ AT₁ Angiotensin II Membrane Preparation

CATALOG NUMBER: HTS064M **QUANTITY:** 200 units
LOT NUMBER: JH1734906 **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: Angiotensin II (Ang II), an octapeptide produced by cleavage of angiotensinogen by angiotensin-converting enzyme, plays a fundamental role in cardiovascular homeostasis. Two GPCRs, AT₁ and AT₂, mediate the effects of AngII, although AT₁ is primarily responsible for the effects of Ang II on renal function and development, thirst, electrolyte and water balance, and arterial blood pressure. Binding of Ang II to AT₁ activates both Gq and Gi (De Gasparo *et al.*, 2000). AT₁ 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 AT₁ interactions and its ligands.

APPLICATIONS: Radioligand Binding Assay

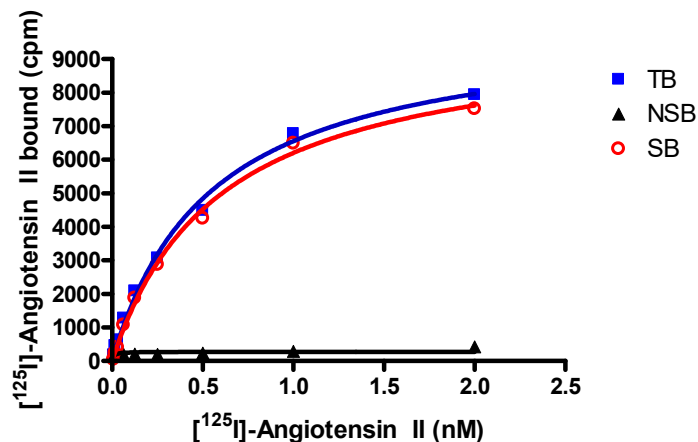


Figure 1. Saturation Binding for AT₁. 5 µg/well AT₁ Membrane Preparation was incubated with increasing amount of [¹²⁵I] labeled Sar¹-Ile⁸-Angiotensin II in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled Angiotensin II. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

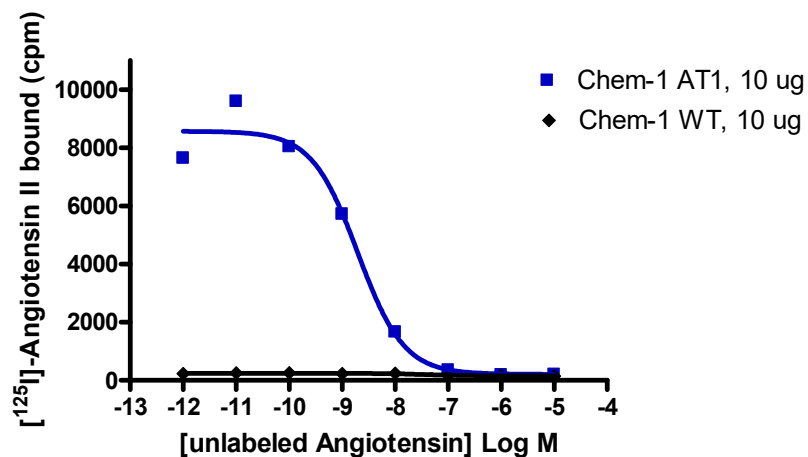


Figure 2. Competition binding for AT₁. AT₁ Membrane Preparation (10 µg/well in a 96-well plate) were incubated with 0.3 nM ¹²⁵I labeled Sar¹-Ile⁸-Angiotensin II and increasing concentrations of unlabeled Angiotensin II, and subjected to filtration binding. Sample data from a representative lot.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} for [¹²⁵I] Angiotensin II binding: 0.45 pmol/mg protein
 K_d for [¹²⁵I] Angiotensin II binding: 0.60 nM

TRANSFECTION: Full-length human AGTR1 cDNA encoding AT₁ (Accession Number: NM_000685).

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous AT₁ 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: [¹²⁵I]-Sar¹-Ile⁸-Angiotensin II. (Perkin Elmer#:NEX-248)

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 40-fold signal:background with ¹²⁵I labeled Sar¹-Ile⁸-Angiotensin II at 0.3 nM.

PRESENTATION: Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA 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. De Gasparo M *et al.* (2000) International Union of Pharmacology. XXIII. The angiotensin II receptors. *Pharmacol. Rev.* 52: 414-472.

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.