

PRODUCT DATASHEET
ChemiScreen™ NK₁ TACHYKININ Membrane Preparation

CATALOG NUMBER:	HTS080M	QUANTITY:	200 units
LOT NUMBER:	21E1807	VOLUME/CONCENTRATION:	1 mL, 2.0 mg/mL

BACKGROUND: The tachykinin peptide family in mammals comprises three peptides, substance P, neurokinin A and neurokinin B, which bind to the tachykinin receptor family of GPCRs, NK₁, NK₂ and NK₃ (Severini *et al.*, 2002). Tachykinins have prominent activity in the GI system, in which they stimulate intestinal contraction and salivation. These effects are mediated by NK₁ and NK₂, and an antagonist of NK₁, aprepitant, is used for treatment of chemotherapy-induced emesis (Rupniak and Kramer, 1999). The NK₁ tachykinin receptor is expressed in brain, and is thought to be involved in depression and nociception (Saria, 1999). Although NK₁ appears to promote nociception in animal models, the significance of this activity is controversial, as antagonists of NK₁ have not proven efficacious in pain relief in humans (Hill, 2000). EMD Millipore's NK₁ 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 NK₁ interactions with Substance P. The membrane preparations exhibit a K_d of 0.28-0.4 nM for [¹²⁵I]-Substance P. With 10 μg/well NK₁ Membrane Prep and 0.25 nM [¹²⁵I]-Substance P, a greater than 5-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand Binding Assay

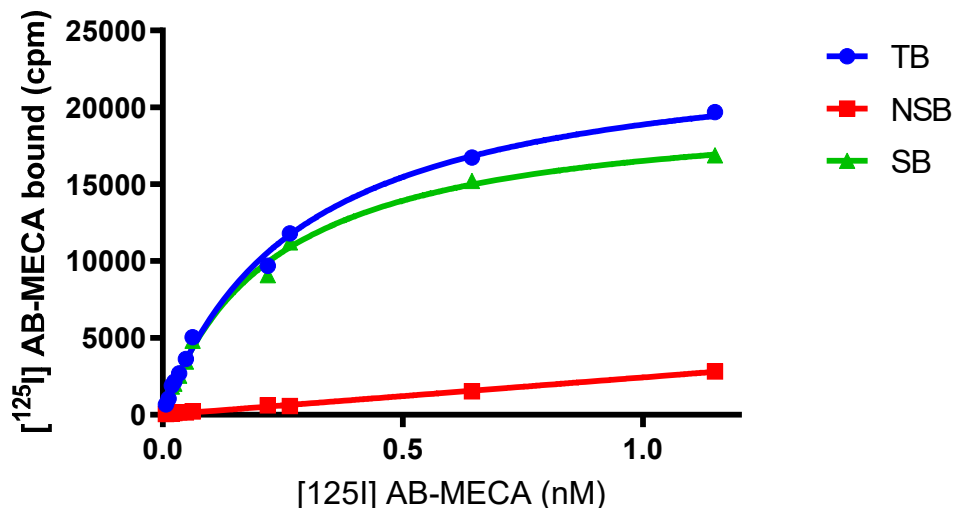


Figure 1. Saturation binding for NK₁. 10 μg/well NK₁ Membrane Preparation was incubated with increasing amount of [¹²⁵I]-Substance P in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 1000-fold excess unlabeled 1 μM [Sar⁹,Met(O₂)¹¹]-substance P.. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

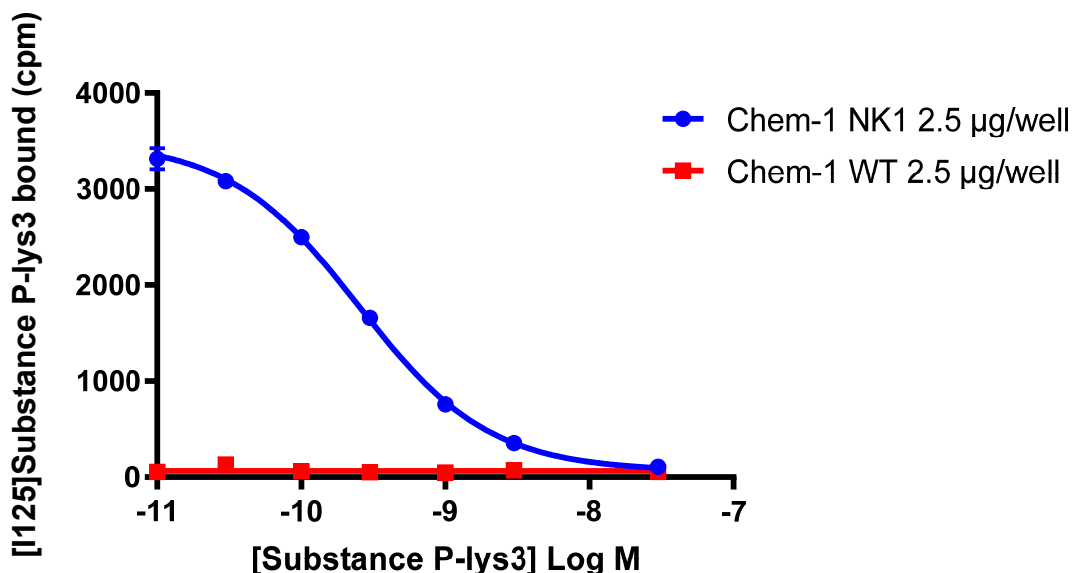


Figure 2. Competition binding for NK₁. NK₁ Membrane Preparation (10 µg/well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 0.25 nM [¹²⁵I]-Substance P and increasing concentrations of unlabeled [Sar⁹,Met(O₂)¹¹]-substance P, and more than 5-fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 10 µg
 B_{max} for [³H] SCH-23390 binding: 1.8 pmol/mg protein
 K_d for [³H] SCH-23390 binding: 0.13 nM
 Signal:background: ≥5-fold

Species: Full-length human TACR1 cDNA encoding NK₁ (Accession Number: NM_001058)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous D₁ 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 50 mM 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 Tris-HCl (pH 7.4), 5 mM MnCl₂, 0.2 % BSA and 40 µg/ml bacitracin, filtered and stored at 4°C.

Radioligand: [¹²⁵I] Substance P (Perkin Elmer # NEX190)

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 10-fold signal:background with [¹²⁵I] Substance P at 0.05 nM.

PRESENTATION:

Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA 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. Avoid repeated freeze/thaw cycles.

REFERENCES:

1. Missale C *et al.* (1998). Dopamine receptors: from structure to function. *Physiol. Rev.* 78:189-225.

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.