

PRODUCT DATASHEET
ChemiScreen™ sst₄ Somatostatin Membrane Preparation

CATALOG NUMBER: HTS125M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: Somatostatin is a 14 or 28 amino acid regulatory peptide that inhibits hormone secretion from the pituitary, pancreas, and other endocrine sites. A family of 6 GPCRs, sst₁, sst_{2A}, sst_{2B}, sst₃, sst₄ and sst₅, mediate the biological activity of somatostatins. The somatostatin receptors couple to G_i to inhibit cAMP production, and also increase MAP kinase signalling. Several tumors have been shown to overexpress somatostatin receptors, and binding of somatostatin to these tumor cells stimulates or inhibits proliferation, depending on the receptor subtypes expressed (Olias *et al.*, 2004). Somatostatin has been implicated in seizure susceptibility in animal models, and activation of sst₄ with selective agonists increases seizure activity (Moneta *et al.*, 2002). sst₄ 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 sst₄ interactions with somatostatin.

APPLICATIONS: Radioligand binding assay

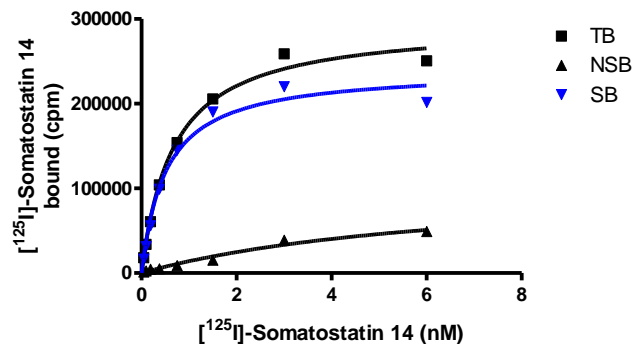


Figure 1. Saturation binding for sst₄. 5 µg/well sst₄ Membrane Preparation was incubated with increasing amount of [¹²⁵I]-Somatostatin 14 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of more than 5000-fold excess unlabeled somatostatin. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

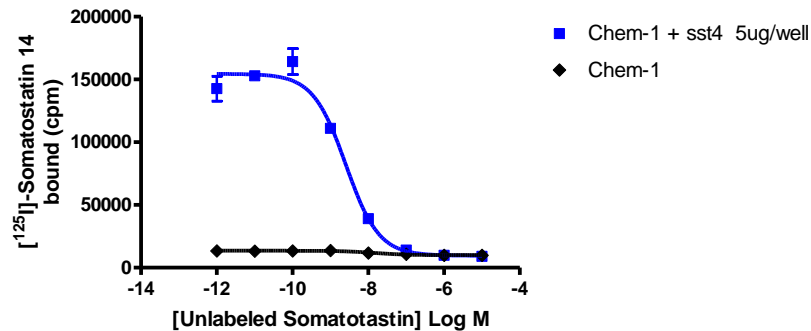


Figure 2. Competition binding for sst₄. sst₄ Membrane Preparation (5 µg/well) or Wild-Type Chem-1 membrane preparation (WT; Catalog # HTS000MC1) was incubated with 1 nM [¹²⁵I]-Somatostatin 14 and increasing concentrations of unlabeled somatostatin, and more than 15-fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 5 µg membrane preparation

B_{max} 20 pmol/mg

K_d 0.5 nM

Signal:background: ≥15-fold

Species: Human SSTR4 encoding SST₄ (Accession Number: NM_001052)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous sst₄ 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]-Somatostatin 14 (Perkin Elmer # NEX389)

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 15-fold signal:background with ¹²⁵I-labeled somatostatin 14 at 1 nM.

PRESENTATION:

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

Packaging method: Membranes protein 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. Moneta D *et al.* (2002) Somatostatin receptor subtypes 2 and 4 affect seizure susceptibility and hippocampal excitatory neurotransmission in mice. *Eur. J. Neurosci.* 16: 843-9.
2. Olias G *et al.* (2004) Regulation and function of somatostatin receptors. *J. Neurochem.* 89: 1057-1091.

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