

PRODUCT DATASHEET

ChemiScreen™ SST₂ Receptor Membrane Preparation

CATALOG NUMBER:	HTS028M	QUANTITY:	200 units
LOT NUMBER:	SC060773	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL

BACKGROUND:

Somatostatin (SST) is a multifunctional peptide with two biologically active forms, SST-14 and SST-28, which are synthesized in neurons throughout the brain as well as in peripheral tissues such as the pancreas and the gut (Gillies, 1997). SST exerts a diverse array of effects that include inhibition of endocrine secretion, modulation of neurotransmission, and regulation of cell proliferation by stimulating a family of five G-protein-coupled receptors. Somatostatin receptor sst₂ mRNA is predominantly expressed in central nervous system. A study using sst₂ knock-out mice has found increased anxiety-related behavior, while locomotor and exploratory activity was decreased in stress-inducing situations (coupled with an increase in pituitary ACTH release, a regulator of the stress response) (Viollet *et al.*, 2000). In the periphery, inhibition of glucagon release by SST in mouse islets is primarily mediated via SST₂ (Strowski *et al.*, 2000). In addition, endogenous SST functions through SST₂ to suppress gastric acid secretion through inhibition of gastrin activity (Martinez *et al.*, 1998). SST₂ 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 of SST₂. The membrane preparations exhibit a K_d of 1.5 nM for [¹²⁵I]-Somatostatin. With 0.75 nM [¹²⁵I]-Somatostatin, 5 µg/well of SST₂ Membrane Prep typically yields greater than a 14-fold signal-to-background ratio.

APPLICATIONS:

Radioligand Binding Assay

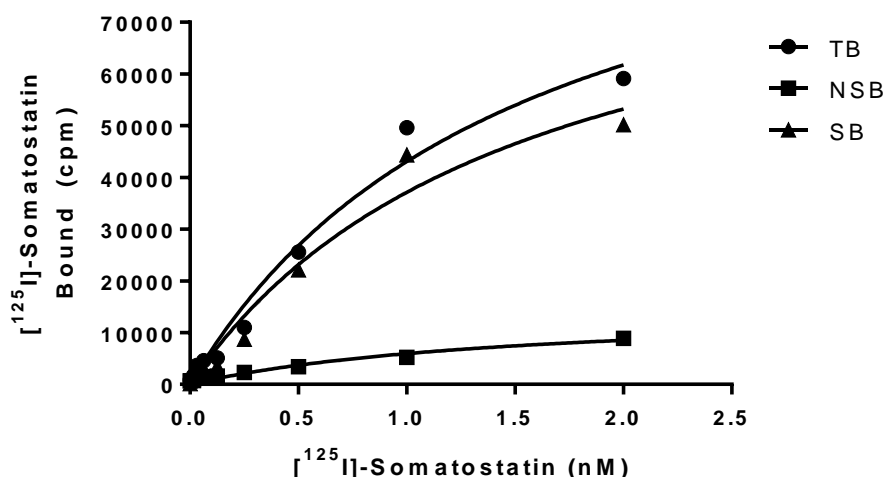


Figure 1. Saturation Binding for SST₂. 5 µg/well of SST₂ Membrane Preparation was incubated with increasing amounts of [¹²⁵I]-Somatostatin in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled somatostatin 14. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample of lot SC060773.

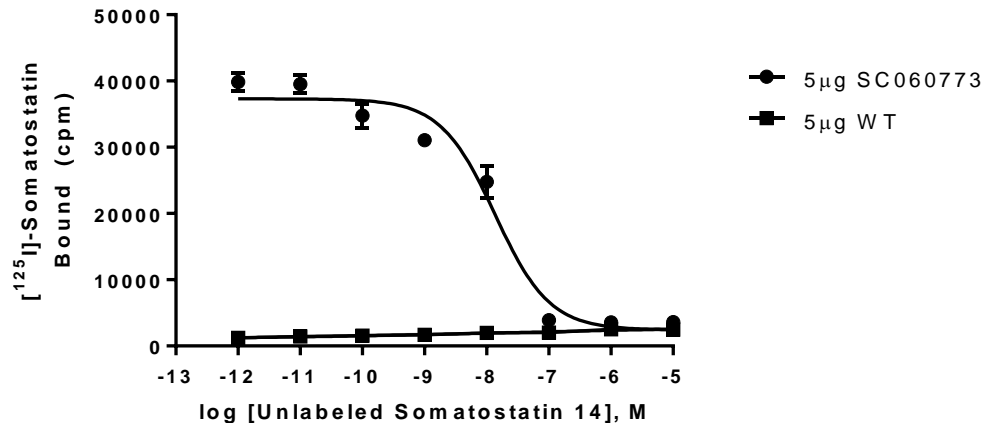


Figure 2. Competition Binding for SST₂. 5 µg/well of SST₂ Membrane Preparation or Wild-type Chem-1 Membrane Preparation (catalog # HTS000MC1) were incubated in a 96-well plate with 0.75 nM [¹²⁵I]-Somatostatin and increasing concentrations of unlabeled somatostatin 14. More than a 14-fold signal:background ratio was obtained. The data are from a representative sample of lot SC060773.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} for [¹²⁵I]-Somatostatin Binding: 7.7 pmol/mg protein
 K_d for [¹²⁵I]-Somatostatin Binding: 1.5 nM
 Signal:background: ≥14-fold

TRANSFECTION: Full-length human sst₂ cDNA encoding sst₂ (Accession Number: NM_001050.2)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous SST₂ 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% polyethylenimine 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 to determine the amount of receptor-associated radioligand binding.

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 (PerkinElmer#: 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 a 14-fold signal:background ratio with [¹²⁵I]-Somatostatin at 0.75 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 1 mg/mL 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. Gillies G (1997). Somatostatin: the neuroendocrine story. *Trends Pharmacol. Sci.* 18:87-95.
 2. Martinez V *et al.* (1998). High basal gastric acid secretion in somatostatin receptor subtype 2 knockout mice. *Gastroenterology* 114:1125-1132.
 3. Strowski MZ *et al.* (2000). Somatostatin inhibits insulin and glucagon secretion via two receptors subtypes: an in vitro study of pancreatic islets from somatostatin receptor 2 knockout mice. *Endocrinology* 141:111-117.
 4. Viollet C *et al.* (2000). Involvement of sst_2 somatostatin receptor in locomotor, exploratory activity and emotional reactivity in mice *Eur. J. Neurosci.* 12:3761-3770.

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.