

PRODUCT DATASHEET
ChemiScreen™ 5-HT₆ Serotonin Membrane Preparation

CATALOG NUMBER: HTS111M **QUANTITY:** 200 units
LOT NUMBER: 2324477 **VOLUME/CONCENTRATION:** 1 mL, 1 mg/mL

BACKGROUND: The neurotransmitter serotonin/5-hydroxytryptamine (5-HT) regulates a wide variety of neurological functions. A family of 13 receptors (12 GPCRs and one ion channel) mediate the effects of serotonin. The serotonin receptor 5-HT₆ is a Gs coupled receptor expressed solely in the CNS, primarily in the limbic and cortical regions. 5-HT₆ appears to play a role in memory and learning, obesity, psychosis, anxiety and epilepsy (Woolley et al., 2004; Fisas et al., 2006). In particular, a 5-HT₆-selective agonist caused significant weight loss in a rat model of diet-induced obesity. 5-HT₆ 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 5-HT₆. The membrane preparations exhibit a K_d of 0.44 nM for [¹²⁵I]-SB258585. With 0.25nM [¹²⁵I]-SB258585, 5µg/well 5-HT₆ Membrane Prep typically yields greater than 6-fold signal-to-background ratio.

APPLICATIONS: Radioligand binding assay

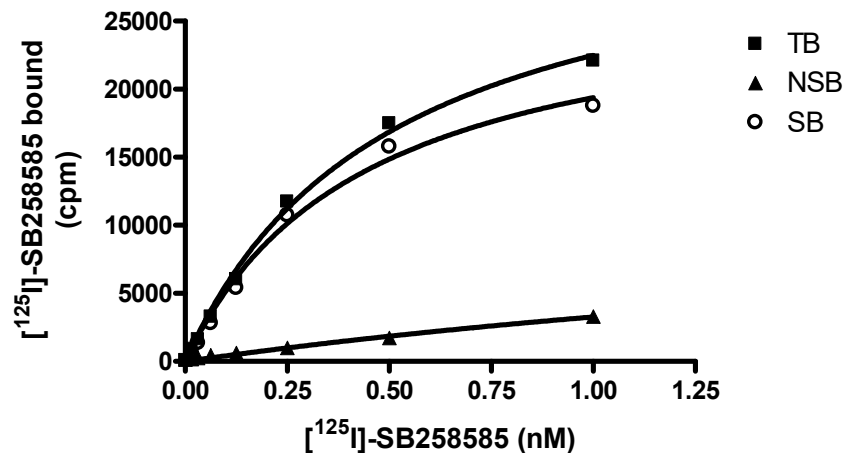


Figure 1. Saturation binding for 5-HT₆. 5 µg/well 5-HT₆ Membrane Preparation was incubated with increasing amount of ¹²⁵I-labeled SB-258585 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled 5-HT. Specific binding (SB) was determined by subtracting NSB from TB.

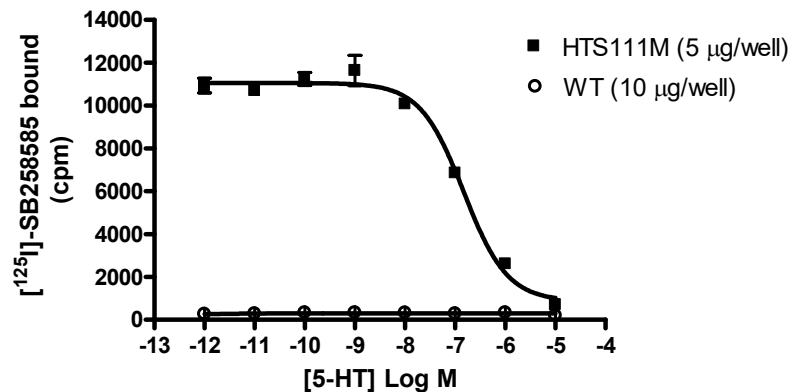


Figure 2. Competition binding for 5-HT₆. 5-HT₆ Membrane Preparation (5 µg/well) and wild-type Chem-1 Membrane Preparation (10 µg/well; Catalog # HTS000MC1) were incubated in a 96-well plate with 0.25 nM ¹²⁵I-labeled SB258585 and increasing concentrations of unlabeled 5-HT. More than 6-fold signal:background was obtained.

SPECIFICATIONS: 1 unit = 5 µg membrane preparation
 B_{max}: 6.1 pmol/mg protein
 K_d: 0.44 nM
 Signal:background: >6

TRANSFECTION: Full-length human HTR6 cDNA encoding the 5-HT₆ Serotonin Receptor (Accession Number: NM_000871)

Species: Human

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous 5-HT₆ 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, then washed with the binding buffer. 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.

Assay Buffer: 20 mM HEPES, 3 mM MgCl₂, 2 mM Ascorbic acid, pH 7.4, filtered and stored at 4°C

Radioligand: [¹²⁵I]-SB258585 (Perkin Elmer #:NEX424)

Wash Buffer: same as binding buffer

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 ¹²⁵I labeled SB258585 at 0.25 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 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. Do not freeze and thaw.

- REFERENCES:**
1. Woolley ML *et al.* (2004) 5-HT₆ receptors. *Curr. Drug Targets CNS Neurol. Disord.* 3: 59-79.
 2. Fisas A *et al.* (2006) Chronic 5-HT₆ receptor modulation by E-6837 induces hypophagia and sustained weight loss in diet-induced obese rats. *Br. J. Pharmacol.* 148: 973-83.

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