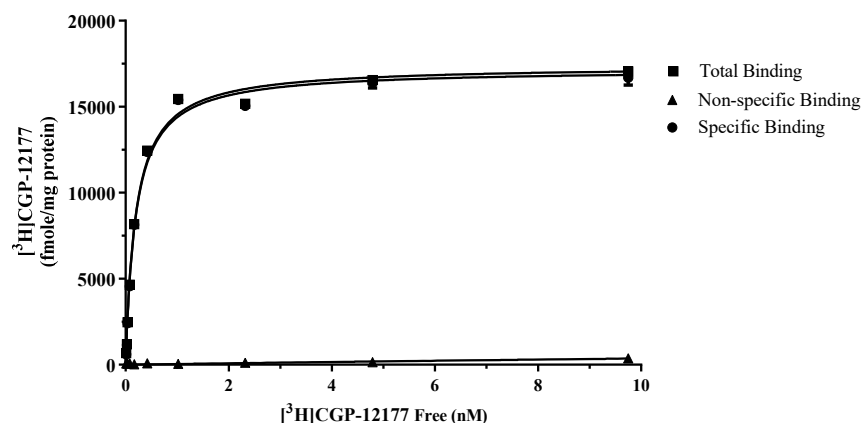


**PRODUCT DATASHEET**
**ChemiScreen™  $\beta_2$  Adrenoceptor Membrane Preparation**

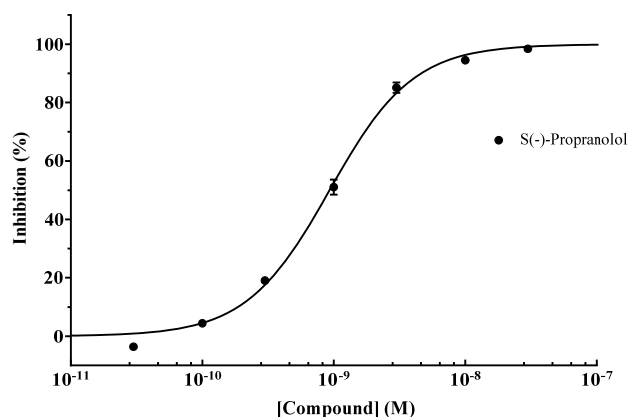
<b>CATALOG NUMBER:</b>	HTS073M	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	22G0603	<b>VOLUME/CONCENTRATION:</b>	1 mL, 1 mg/mL

**BACKGROUND:** The endogenous catecholamines epinephrine and norepinephrine have profound effects on smooth muscle activity, cardiac function, carbohydrate and fat metabolism, hormone secretion, neurotransmitter release, and central nervous system actions. These activities are mediated by GPCRs belonging to two subfamilies, the  $\alpha$ - and  $\beta$ -adrenergic receptors (Bylund *et al.*, 1994). The  $\beta$ -adrenergic receptors, primarily the  $\beta_2$  subtype, mediate relaxation of smooth muscle in many tissues, and  $\beta_2$ -selective agonists are the preferred drugs for stimulating bronchodilation in the treatment of asthma and chronic obstructive pulmonary disease (Sears and Lotvall, 2005). Activation of the  $\beta$ -adrenergic receptors, primarily the  $\beta_1$  subtype and to a lesser extent the  $\beta_2$  subtype, acutely increases heart rate, cardiac output, and cardiac automaticity, and chronically increases cardiac myocyte apoptosis. As a result,  $\beta$ -adrenergic receptor antagonists ( $\beta$  blockers) are effective in the treatment of congestive heart failure and arrhythmia (Feldman *et al.*, 2005). The  $\beta_2$  adrenoceptor 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  $\beta_2$  adrenoceptor interactions with (-)-ludocyanopindolol (ICYP). The membrane preparations exhibit  $K_d$  of 0.195 nM for [ $^3$ H]CGP-12177. With 5  $\mu$ g/well of  $\beta_2$  Adrenoceptor Membrane Prep and 0.2 nM [ $^3$ H]CGP-12177, a greater than 10-fold signal-to-background ratio was obtained.

**APPLICATIONS:** Radioligand Binding Assay



**Figure 1. Saturation Binding for  $\beta_2$  Adrenoceptor.** 5  $\mu$ g/well of  $\beta_2$  Adrenoceptor Membrane Preparation was incubated with increasing amounts of [ $^3$ H]CGP-12177 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled ICI 118,551. Specific binding (SB) was determined by subtracting NSB from TB.



**Figure 2. Competition binding for  $\beta_2$  adrenoceptor.** 5  $\mu\text{g}$ /well of  $\beta_2$  adrenoceptor Membrane Preparation were incubated with 0.2 nM [ $^3\text{H}$ ]CGP-12177 and increasing concentrations of (S)-(-)-Propranolol, and more than a 10-fold signal:background was obtained.

**SPECIFICATIONS:** 1 unit = 5  $\mu\text{g}$  membrane preparation  
 B<sub>max</sub> for [ $^3\text{H}$ ]CGP-12177 Binding: 17.2 pmol/mg  
 K<sub>d</sub> for [ $^3\text{H}$ ]CGP-12177 Binding: 0.195 nM  
 Signal:Background: >10-fold

**Species:** Full length human ADRB2 encoding  $\beta_2$  adrenoceptor (Accession number NM\_000024).

**HOST CELLS:** Chem-2, a suspension mammalian cell line without any endogenous  $\beta_2$  adrenoceptor 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% polyethyleneimine for 30 min, then washed with 50 mM HEPES, pH 7.4. 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 MgCl<sub>2</sub>, 1 mM CaCl<sub>2</sub>, filtered and stored at 4°C

**Radioligand:** [ $^3\text{H}$ ] CGP-12177 (PerkinElmer, NET-1061)

**Wash Buffer:** 50 mM HEPES, pH 7.4, 500 mM NaCl, 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 40-fold signal:background ratio with [ $^{125}\text{I}$ ]-(-)Iodocyanopindolol at 0.5 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 in 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. Bylund DB *et al.* (1994). IV. International Union of Pharmacology nomenclature of adrenoceptors. *Pharmacol. Rev.* 46:121-136.
2. Feldman DS *et al.* (2005). Mechanisms of Disease:  $\beta$ -adrenergic receptors—alterations in signal transduction and pharmacogenomics in heart failure. *Nat. Clin. Pract. Cardiovasc. Med.* 2:475-83.
3. Sears MR and Lotvall J (2005). Past, present and future— $\beta_2$ -adrenoceptor agonists in asthma management. *Respir. Med.* 99:152-70.

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.