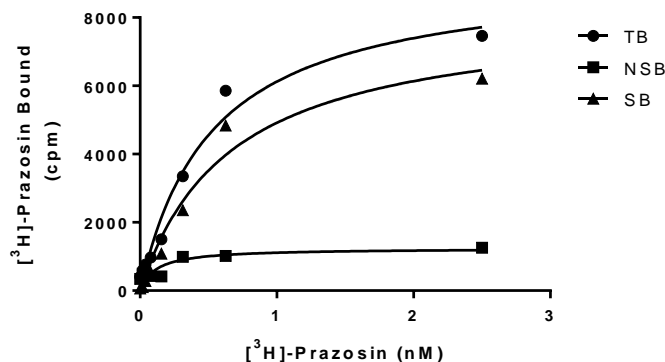


**PRODUCT DATASHEET**
**ChemiScreen™  $\alpha_{1B}$  Adrenergic Membrane Preparation**

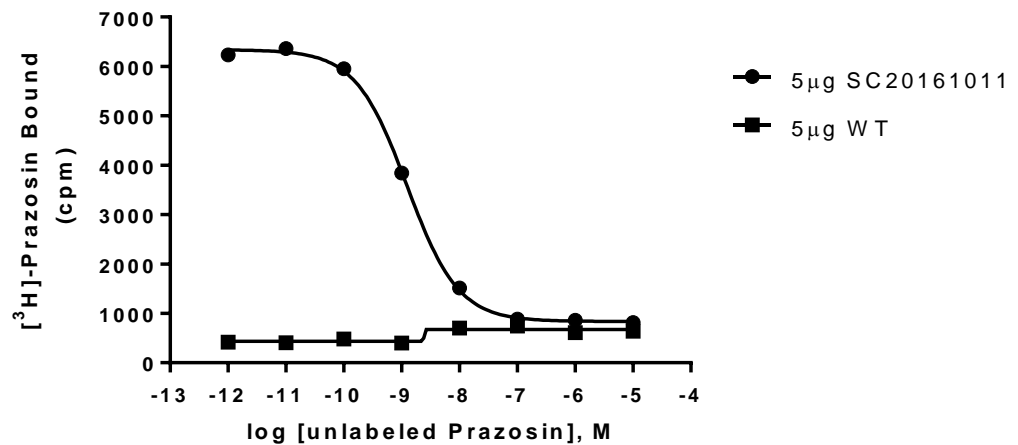
<b>CATALOG NUMBER:</b>	HTS158M	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	SC20161011	<b>VOLUME/CONCENTRATION:</b>	1 mL, 1 mg/mL

**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  $\beta$ -adrenoceptors (Bylund *et al.*, 1994). The three members of the  $\alpha_1$  subclass of adrenoceptors,  $\alpha_{1A}$ ,  $\alpha_{1B}$  and  $\alpha_{1D}$ , couple to  $G_q$ , and promote contraction of vascular and urinary tract smooth muscle, relaxation of intestinal smooth muscle, increased contractile force in the heart, and glycogenolysis and gluconeogenesis in the liver. The different subtypes have overlapping distributions and variably contribute to these effects depending on species and tissue. Overexpression of a constitutively active  $\alpha_{1B}$  mutant in the heart of transgenic mice resulted in cardiac hypertrophy with increased heart weight/body weight ratios. Analysis of  $\alpha_{1B}$  knock out mice has provided evidence that  $\alpha_{1B}$  is a mediator of blood pressure and aortic contractile responses induced by  $\alpha_1$  agonists (Milano *et al.*, 1994). The locomotor and rewarding effects of psychostimulants and opiates were suppressed in mice lacking  $\alpha_{1B}$ -adrenergic receptors (Drouin *et al.* 2002).  $\alpha_{1B}$  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  $\alpha_{1B}$ . The membrane preparations exhibit a  $K_d$  of 0.62 nM for [ $^3$ H]-Prazosin. With 1 nM [ $^3$ H]-Prazosin, 5  $\mu$ g/well of  $\alpha_{1B}$  Membrane Prep yields greater than a 5-fold signal-to-background ratio.

**APPLICATIONS:** Radioligand Binding Assay



**Figure 1. Saturation Binding for  $\alpha_{1B}$ .** 5  $\mu$ g/well  $\alpha_{1B}$  Membrane Preparation was incubated with increasing amounts of [ $^3$ H]-Prazosin in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 200-fold excess unlabeled Prazosin. Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample lot.



**Figure 2. Competition Binding for  $\alpha_{1B}$ .** 5 $\mu$ g/well  $\alpha_{1B}$  Membrane Preparation and Wild-type Chem-1 Membrane Preparation (catalog # HTS000MC1) were incubated in a 96-well plate with 1 nM [<sup>3</sup>H]-Prazosin and increasing concentrations of unlabeled Prazosin. More than a 5-fold signal:background ratio was obtained. The data are from a representative sample lot.

**SPECIFICATIONS:** 1 unit = 5  $\mu$ g  
 $B_{max}$  for [<sup>3</sup>H]-Prazosin Binding: 15.1 pmol/mg protein  
 $K_d$  for [<sup>3</sup>H]-Prazosin Binding: 0.62 nM  
Signal:background: >5-fold

**TRANSFECTION:** Full-length human ADRA1B cDNA encoding  $\alpha_{1B}$  adrenergic receptor (Accession Number: NM\_000679.3)

**HOST CELLS:** Chem-1, an adherent mammalian cell line without any endogenous  $\alpha_{1B}$  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 non-binding 96-well plate, and incubated for 2 h. Prior to filtration, an FC 96-well harvest plate is coated with 0.33% polyethyleneimine for 30 min, and washed with 50 mM Tris, pH 7.4. The binding reactions are transferred to the filter plate, and washed 3 times (1 mL per well per wash) with Wash Buffer. The wells are then dried and counted.

**Binding Buffer:** 50 mM Tris, pH 7.4, 10 mM MgCl<sub>2</sub>, 1 mM EDTA, filtered and stored at 4°C

**Radioligand:** [<sup>3</sup>H]-Prazosin (PerkinElmer # NET823)

**Wash Buffer:** 50 mM Tris, pH 7.4, 500 mM 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 5-fold signal:background with 1 nM [<sup>3</sup>H]-Prazosin.

**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 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. Bylund DB *et al.* (1994). IV. International Union of Pharmacology nomenclature of adrenoceptors. *Pharmacol. Rev.* 46:121-136.
2. Cavalli A *et al.* (1997). Decreased blood pressure response in mice deficient of the  $\alpha_{1B}$ -AR. *Proc. Natl. Acad. Sci. USA* 94:11589–11594.
3. Milano CA *et al.* (1994). Myocardial expression of a constitutively active  $\alpha_{1B}$ -adrenergic receptor in transgenic mice induces cardiac hypertrophy. *Proc. Natl. Acad. Sci. USA* 91:10109-10113.

**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.