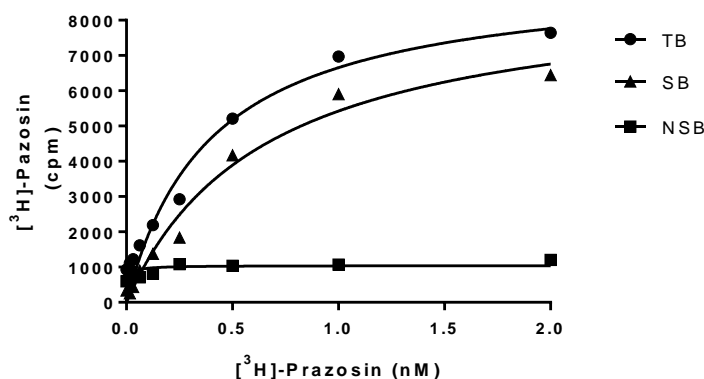


**PRODUCT DATASHEET**
**ChemiScreen™  $\alpha_{1D}$  Adrenergic Membrane Preparation**

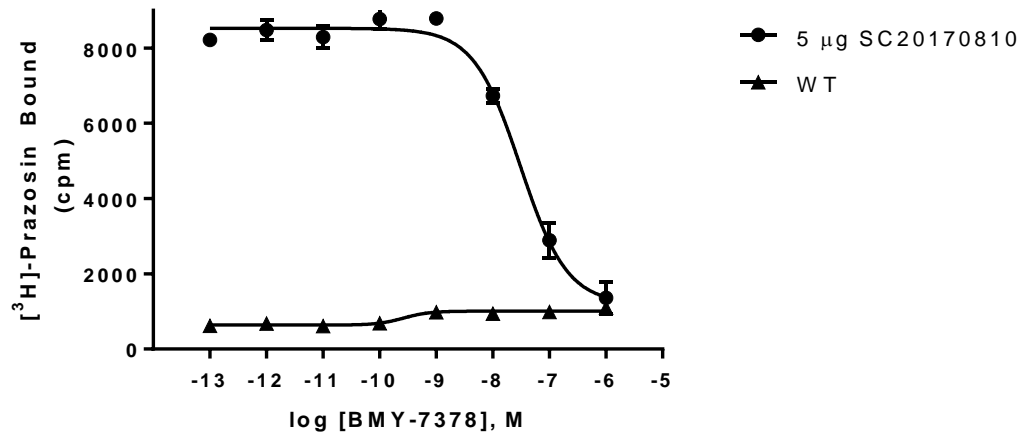
<b>CATALOG NUMBER:</b>	HTS216M	<b>QUANTITY:</b>	200 units
<b>LOT NUMBER:</b>	SC20170810	<b>VOLUME/CONCENTRATION:</b>	1 mL, 1 mg/mL

**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. The  $\alpha_{1D}$  adrenergic receptor mediates smooth muscle contraction in several tissues. In the vasculature, activation of  $\alpha_{1D}$  increases blood pressure (Tanoue *et al.*, 2002; Hosoda *et al.*, 2005). In the urinary tract,  $\alpha_{1D}$  promotes bladder contraction. Antagonists of  $\alpha_1$  receptors are used to treat bladder outlet obstruction, and this effect is thought to be mediated by  $\alpha_{1D}$  (Chen *et al.*, 2005). The  $\alpha_{1D}$  adrenergic receptors has a relatively long N-terminal extracellular domain, and truncation of this domain has been shown to increase expression of the receptor at the cell surface (Pupo *et al.*, 2003). The  $\alpha_{1D}$  membrane preparations, which contain a version of  $\alpha_{1D}$  lacking residues 2-79, 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  $\alpha_{1D}$ .

**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 [<sup>3</sup>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  $\alpha_{1D}$ .**  $\alpha_{1D}(\Delta^{2-79})$  Membrane Preparation (5 and 10  $\mu\text{g/well}$ ) and wild-type Chem-1 Membrane Preparation (10  $\mu\text{g/well}$ , catalog # HTS000MC1) were incubated in a 96-well plate with 0.5 nM [<sup>3</sup>H]-labeled prazosin and increasing concentrations of unlabeled BM Y-7378. More than 5-fold signal:background was obtained

**SPECIFICATIONS:** 1 unit = 5  $\mu\text{g}$   
 $B_{\text{max}}$ : 19.17 pmol/mg  
 $K_{\text{d}}$ : ~0.65 nM  
Signal:background:  $\geq 5$ -fold

**TRANSFECTION:** Truncated human ADRA1D cDNA encoding  $\alpha_{1D}$  lacking residues 2-79 (based on Accession Number: NM\_000678)

**HOST CELLS:** Chem-1, an adherent mammalian cell line without any endogenous  $\alpha_{1D}$  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 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  $\text{MgCl}_2$ , 1 mM  $\text{CaCl}_2$ , 0.2% BSA, filtered and stored at 4°C

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

**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 5-fold signal:background with <sup>3</sup>H labeled prazosin 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 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:**

Bylund DB *et al.* (1994) IV. International Union of Pharmacology nomenclature of adrenoceptors. *Pharmacol. Rev.* 46: 121-136.

Chen Q *et al.* (2005) Function of the lower urinary tract in mice lacking  $\alpha_{1D}$ -adrenoceptor. *J. Urol.* 174: 370-374.

Horie K *et al.* (1995) Selectivity of the imidazoline  $\alpha$ -adrenoceptor agonists (oxymetazoline and cirazoline) for human cloned  $\alpha_1$ -adrenoceptor subtypes. *Br. J. Pharmacol.* 116: 1611-1618.

Hosoda C *et al.* (2005) Two  $\alpha_1$ -adrenergic receptor subtypes regulating the vasopressor response have differential roles in blood pressure regulation. *Mol. Pharmacol.* 67: 912-922.

Pupo AS *et al.* (2003) N-terminal truncation of human  $\alpha_{1D}$ -adrenoceptors increases expression of binding sites but not protein. *Eur. J. Pharmacol.* 462: 1-8.

Tanoue A *et al.* (2002) The  $\alpha_{1D}$ -adrenergic receptor directly regulates arterial blood pressure via vasoconstriction. *J. Clin. Invest.* 109: 765-775.

**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.