

PRODUCT DATASHEET
ChemiScreen™ VPAC₂ VIP/PACAP Membrane Preparation

CATALOG NUMBER:	HTS079M	QUANTITY:	200 units
LOT NUMBER:	SC191274	VOLUME/CONCENTRATION:	1 mL, 1 mg/mL

BACKGROUND: Vasoactive intestinal peptide (VIP), a 28 amino acid peptide originally isolated by its vasodilation activity, binds to two class B GPCRs, VPAC₁ and VPAC₂, to exert its functions in the CNS, vasculature, immune system and adrenal medulla (Harmar *et al.*, 1998). In the immune system, VPAC₂ is expressed on stimulated CD4 T cells, and binding of T cell-derived VIP to VPAC₂ induces a shift toward the Th2 pathway (Delgado *et al.*, 2004; Voice *et al.*, 2004). In addition, VPAC₂ is an essential regulator of the circadian pacemaker of the hypothalamic suprachiasmatic nuclei (Hughes *et al.*, 2004). VPAC₂ 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 VPAC₂ interactions with PACAP. The VPAC₂ receptor membrane preparations exhibit a K_d of 0.86 nM for [¹²⁵I]-PACAP27. With 0.5 nM [¹²⁵I]-PACAP27, 5 µg/well of VPAC₂ Receptor Membrane Prep yields greater than a 10-fold signal-to-background ratio.

APPLICATIONS: Radioligand Binding Assay

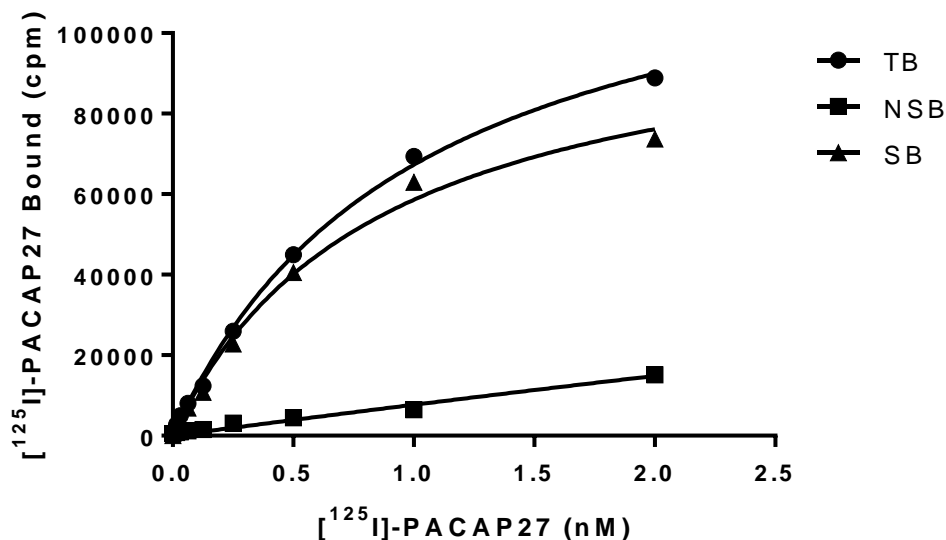


Figure 1. Saturation Binding for VPAC₂. 5 µg/well of VPAC₂ Membrane Preparation were incubated with increasing amounts of [¹²⁵I]-PACAP27 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of 500-fold excess unlabeled PACAP(1-27). Specific binding (SB) was determined by subtracting NSB from TB. The data are from a representative sample of lot SC191274.

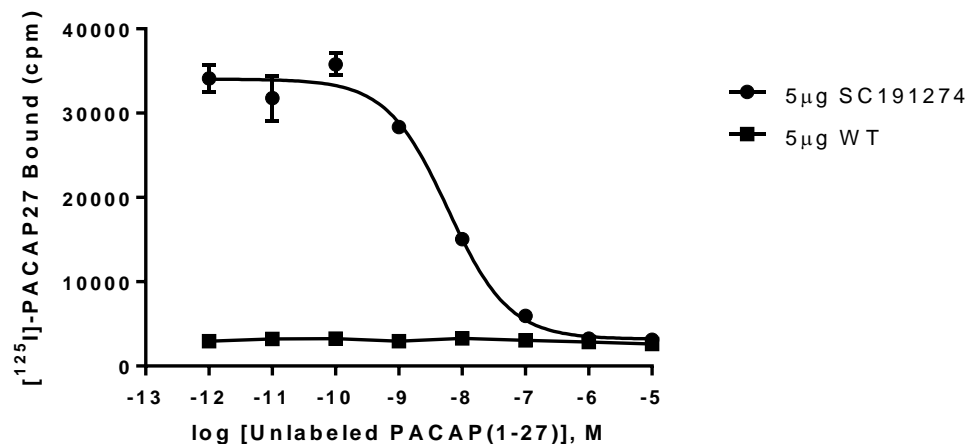


Figure 2. Competition Binding for VPAC₂. 5 µg/well of VPAC₂ Membrane Preparation or Wild-Type Chem-1 membrane preparation (Catalog # HTS000MC1) were incubated with 0.5 nM [¹²⁵I]-PACAP27 and increasing concentrations of unlabeled PACAP(1-27), and more than a 10-fold signal:background ratio was obtained. The data are from a representative sample of lot SC191274.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} for [¹²⁵I]-PACAP27 Binding: 8.9 pmol/mg protein
 K_d for [¹²⁵I]-PACAP27 Binding: 0.86 nM
 Signal:Background: ≥10-fold

Species: Human VPAC₂ (Accession number NM_003382)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous VPAC₂ 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, 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 for determination 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]-PACAP27 (PerkinElmer # NEX294)

Wash Buffer: 50 mM HEPES, 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 10-fold signal:background ratio with [¹²⁵I]-PACAP27 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. Delgado M *et al.* (2004). The significance of vasoactive intestinal peptide in immunomodulation. *Pharmacol. Rev.* 56:249-290.
 2. Harmar AJ *et al.* (1998). International Union of Pharmacology. XVIII. Nomenclature of receptors for vasoactive intestinal peptide and pituitary adenylate cyclase-activating polypeptide. *Pharmacol. Rev.* 50:265-270.
 3. Hughes AT *et al.* (2004). Aberrant gating of photic input to the suprachiasmatic circadian pacemaker of mice lacking the VPAC₂ receptor. *J. Neurosci.* 24:3522-6.
 4. Voice J *et al.* (2004). c-Maf and JunB mediation of Th2 differentiation induced by the type 2 G Protein-Coupled Receptor (VPAC₂) for vasoactive intestinal peptide. *J. Immunol.* 172:7289-7296.

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