

PRODUCT DATASHEET
ChemiScreen™ V_{1B} Vasopressin Membrane Preparation

CATALOG NUMBER: HTS136M **QUANTITY:** 200 units
LOT NUMBER: **VOLUME/CONCENTRATION:** 1 mL, 2 mg/mL

BACKGROUND: Arginine vasopressin (AVP) is a 9 amino acid peptide that functions as an antidiuretic, vasoconstrictor and neurotransmitter. The three vasopressin receptors, V_{1A}, V_{1B} and V₂, are GPCRs; V_{1A} and V_{1B} couple to G_q and calcium release, whereas V₂ couples to G_s (Birnbaumer, 2000). The V_{1B} receptor is expressed prominently in the anterior pituitary, where it mediates vasopressin-induced release of ACTH (Tanoue *et al.*, 2004). A selective antagonist of V_{1B} has recently been developed and shown to reduce depression, anxiety, and aggression in rodents (Blanchard *et al.*, 2005; Griebel *et al.*, 2002). Cloned human V_{1B}-expressing cell line is made in the Chem-1 host, which supports high levels of recombinant V_{1B} expression on the cell surface and contains high levels of the promiscuous G protein G α 15 to couple the receptor to the calcium signaling pathway. V_{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 antagonists of V_{1B} interactions with vasopressin. The membrane preparations exhibit a K_d of 1.03 nM for [³H]-vasopressin. With 10 μ g/well V_{1B} Membrane Prep and 1.5 nM [³H]-vasopressin, a greater than 4-fold signal-to-background ratio was obtained.

APPLICATIONS: Radioligand binding assay

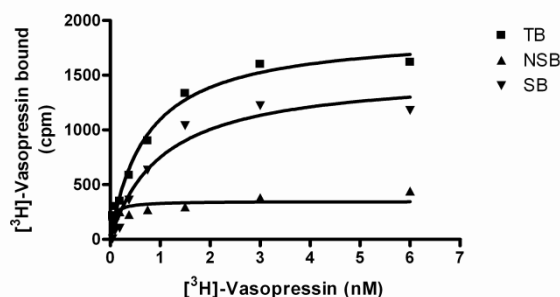


Figure 1. Saturation binding for V_{1B}. 5 μ g/well V_{1B} Membrane Preparation was incubated with increasing amount of [³H]-vasopressin in the absence (total binding, TB) or presence (nonspecific binding, NSB) of greater than 500-fold excess unlabeled vasopressin. Specific binding (SB) was determined by subtracting NSB from TB. Sample data from a representative lot.

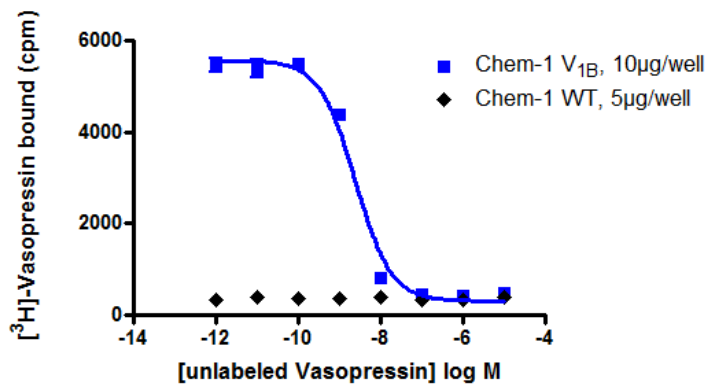


Figure 2. Competition binding for V_{1B}. V_{1B} Membrane Preparation (10 µg/well) or Wild-Type Chem-1 membrane preparation (5 µg/well; Catalog # HTS000MC1) was incubated with 1.5 nM [³H]-vasopressin and increasing concentrations of unlabeled vasopressin, and more than 4-fold signal:background was obtained. Representative sample data.

SPECIFICATIONS: 1 unit = 10 µg membrane preparation
 B_{max}: 2.79 pmol/mg
 K_d: 1.03 nM
 Signal:background: ≥4-fold

Species: Human V_{1B} (Accession number NM_000707)

HOST CELLS: Chem-1, an adherent mammalian cell line without any endogenous V_{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 nonbinding 96-well plate, and incubated for 1-2 h. Prior to filtration, a GF/C 96-well filter plate 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 MgCl₂, 1 mM CaCl₂, 0.2% BSA, filtered and stored at 4°C

Radioligand: [³H] vasopressin (Perkin Elmer # NET800)

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 4-fold signal:background with ³H-labeled vasopressin at 1.5 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 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. Birnbaumer M (2000) Vasopressin receptors. *Trends Endocrinol. Metab.* 11:406-10.
2. Blanchard RJ *et al.* (2005) AVP V1b selective antagonist SSR149415 blocks aggressive behaviors in hamsters. *Pharmacol. Biochem. Behav.* 80: 189-194.
3. Griebel G *et al.* (2002) Anxiolytic- and antidepressant-like effects of the non-peptide vasopressin V_{1b} receptor antagonist, SSR149415, suggest an innovative approach for the treatment of stress-related disorders. *Proc. Natl. Acad. Sci. USA* 99: 6370-6375.
4. Tanoue A *et al.* (2004) The vasopressin V1b receptor critically regulates hypothalamic-pituitary-adrenal axis activity under both stress and resting conditions. *J. Clin. Invest.* 113: 302-309.

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