



# Product Manual

**Modified Procedure!  
Read Entire Manual  
Prior to Using Kit!**

## **Arg<sup>8</sup>-Vasopressin ELISA kit**

Catalog #: ADI-901-017A

5x96 Well Kit



## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Do not mix components from different kit lots or use reagents beyond the expiration date of the kit.



Please contact Enzo Life Sciences Technical Support if necessary.

## TABLE OF CONTENTS

Description .....	2
Introduction .....	2
Principle .....	3
Materials Supplied .....	4
Storage .....	5
Other Materials Needed.....	5
Sample Handling.....	6
Reagent Preparation.....	8
Assay Procedure.....	9
Calculation of Results .....	10
Typical Results.....	11
Typical Standard Curve .....	11
Performance Characteristics.....	12
Sample Dilution Recommendations .....	15
References.....	16
Contact Information.....	18



For proper performance, use the insert provided with each individual kit received.

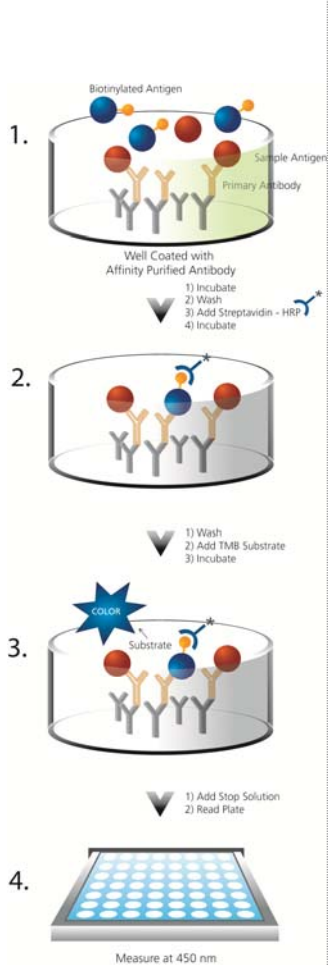
## DESCRIPTION

The Arg<sup>8</sup>-Vasopressin ELISA kit is a competitive immunoassay for the quantitative determination of Vasopressin in plasma, serum, and conditioned cell media. Please read the entire kit manual before performing the assay. For further explanation of the principles and practice of immunoassays please see the excellent books by Chard<sup>1</sup> or Tijssen<sup>2</sup>.

## INTRODUCTION

Arginine Vasopressin (AVP) is a 9 amino acid peptide with a 6-member disulfide ring. It is structurally related to oxytocin differing in 2 amino acids. It is synthesized in the hypothalamus supraoptic and paraventricular nuclei. It is stored in the posterior pituitary for release. AVP has powerful antidiuretic action and is also known as antidiuretic hormone (ADH)<sup>2</sup>. It acts upon the collecting tubule of the kidney increasing permeability to water and urea. It also has neurotransmitter and peripheral humoral functions. AVP has been shown to be released upon both osmotic and non-osmotic stimuli<sup>4,5</sup>, and its release into peripheral blood causes effects upon a number of factors, including emotional stress, posture, blood volume, and temperature<sup>6-9</sup>. Alcohol appears to inhibit AVP secretion. Serum AVP measurement is used clinically for studies involving diabetes insipidus, syndrome of inappropriate ADH secretion (SIADH), ectopic AVP production and psychogenic water intoxication<sup>10-14</sup>.

## PRINCIPLE



1. Standards and samples are added to wells coated with a goat-anti-rabbit IgG antibody. A blue solution of Arg<sup>8</sup>-Vasopressin conjugated to biotin is added, followed by a yellow solution of a rabbit polyclonal antibody to Arg<sup>8</sup>-Vasopressin.
2. The plate is incubated overnight at 4°C. During this incubation the specific antibody binds, in a competitive manner, the Arg<sup>8</sup>-Vasopressin in the sample or conjugate. The plate is then washed, leaving only bound Arg<sup>8</sup>-Vasopressin or the biotinylated Vasopressin conjugate.
3. A solution of streptavidin conjugated to horseradish peroxidase is added to each well, to bind the biotinylated Arg<sup>8</sup>-Vasopressin. The plate is incubated shaking at room temperature.
4. The plate is washed to remove unbound HRP conjugate. TMB substrate solution is added. An HRP-catalyzed reaction generates a blue color in the solution.
5. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at **450nm**. The amount of signal is inversely proportional to the concentration of Arg<sup>8</sup>-Vasopressin in the sample.

## MATERIALS SUPPLIED



The standard should be handled with care due to the known and unknown effects of the antigen.



Activity of conjugate is affected by nucleophiles such as azide, cyanide, and hydroxylamine.



Protect TMB substrate from prolonged exposure to light.



Stop Solution 2 is caustic. Keep tightly capped.

- 1. Goat anti-Rabbit IgG Microtiter Plate, Five Plates of 96 Wells**  
**Catalog No. 80-0060**  
Plates using break-apart strips coated with goat antibody specific to rabbit IgG.
- 2. Assay Buffer 28, 100ml**  
**Catalog No. 80-1797**  
Phosphate buffered saline containing BSA and detergent.
- 3. Arg<sup>8</sup>-Vasopressin Standard, 3 x 0.5ml**  
**Catalog No. 80-2698**  
A solution of 10,000pg/ml Vasopressin.
- 4. Arg<sup>8</sup>-Vasopressin Conjugate, 25ml**  
**Catalog No. 80-2700**  
A blue solution of biotin conjugated with Vasopressin.
- 5. Arg<sup>8</sup>-Vasopressin Antibody, 25ml**  
**Catalog No. 80-2702**  
A yellow solution of a rabbit polyclonal antibody to Vasopressin.
- 6. Wash Buffer Concentrate, 100ml**  
**Catalog No. 80-1287**  
Tris buffered saline containing detergents.
- 7. SA-HRP, 100ml**  
**Catalog No. 80-2704**  
A solution of streptavidin conjugated to horseradish peroxidase.
- 8. TMB Substrate, 100ml**  
**Catalog No. 80-2035**  
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide.
- 9. Stop Solution 2, 50ml**  
**Catalog No. 80-1804**  
A 1N solution of hydrochloric acid in water.
- 10. Vasopressin Assay Layout Sheet, 1 each**  
**Catalog No. 30-0041**
- 11. Plate Sealer, 5 each**  
**Catalog No. 30-0012**



Reagents  
require  
separate  
storage  
conditions.

## STORAGE

All components of this kit, **except the standard**, are stable at 4°C until the kit's expiration date. Upon receipt, the standard must be stored frozen at -20°C. Storage conditions do not necessarily reflect shipping conditions.

## OTHER MATERIALS NEEDED

1. Deionized or distilled water.
2. Precision pipets for volumes between 5µl and 1,000µl.
3. Repeater pipets for dispensing 50µl and 200µl.
4. Disposable beaker for diluting buffer concentrates.
5. Graduated cylinders.
6. Lint-free paper for blotting.
7. Microplate reader capable of reading at 450nm.
8. A microplate shaker.
9. Nitrogen gas (required for the suggested extraction protocol)
10. Acetone (required for the suggested extraction protocol)
11. Petroleum Ether (required for the suggested extraction protocol)
12. Butanol (optional for the suggested extraction protocol)
13. Diisopropyl ether (optional for the suggested extraction protocol)

## SAMPLE HANDLING



Sample handling procedures should be completed prior to reagent preparation.

The Arg<sup>8</sup>-Vasopressin ELISA kit is compatible with Vasopressin samples in a number of matrices. Vasopressin samples should be reconstituted in kit Assay Buffer 28 for extracted serum and plasma samples and diluted into Assay Buffer 28 for tissue culture medium samples. Please refer to the Sample Recovery recommendations on page 15 for details of suggested dilutions. However, the end user must verify that the recommended dilutions are appropriate for their samples. **Samples containing rabbit IgG may interfere with the assay.**

Due to the low endogenous levels of native Arg<sup>8</sup>-Vasopressin we recommend extraction of samples, thereby concentrating them and allowing for accurate determinations of Arg<sup>8</sup>-Vasopressin. An extraction protocol is outlined below. Because of the labile nature of Vasopressin we recommend several precautions in collecting and analyzing samples.



Samples must be stored at or below -20°C to avoid loss of bioactive analyte. Avoid repeated freeze/thaw cycles.

Blood samples should be drawn into chilled EDTA (1mg/ml blood) or serum tubes containing Aprotinin (500 KIU/ml of blood). Centrifuge the samples at 1,600 x g for 15 minutes at 4°C. Transfer the plasma or serum to a plastic tube and store at -70°C or lower for long term storage. Avoid repeated freeze/thaw cycles. The stability of some peptides is improved by the addition of a protease inhibitor cocktail to the sample before freezing.

If samples are thought to be lipemic, the following procedure can be used to delipidate prior to extraction.

1. Prepare mixture of 40:60 butanol:diisopropyl ether. Vortex.
2. Add equal volume of butanol:diisopropyl ether to sample. Vortex.
3. Centrifuge at 8,000 x g for 5 minutes.
4. Remove top organic layer and discard. Measure aqueous layer and transfer to new tube.

### Extraction Procedure:

1. Add 2x volume of ice cold acetone to sample. Vortex.
2. Centrifuge at 3,000 x g for 20 minutes.
3. Transfer supernatant to new tube.
4. Add 5x volume of ice cold petroleum ether. Vortex.
5. Centrifuge at 3,000 x g for 10 minutes.
6. Discard top ether layer. Carefully transfer remaining aqueous layer to glass tube and dry down under gas.
7. Reconstitute sample with Assay Buffer.



If the end user chooses to vary from the extraction procedure noted, it is up to the end user to determine the appropriate dilution of samples and assay validation.



Please note that recovery of peptides from extraction processes can be variable. It is important to optimize any process to obtain optimum recoveries. Extraction efficiencies can be determined by a number of methods, such as spiking into paired samples and determining the recovery of this known amount of added Vasopressin.

Extraction efficiencies for matrices tested are listed below. For each matrix listed, high, middle, and low concentrations of Arg<sup>8</sup>-Vasopressin were spiked into the matrix, then extracted as per the sample extraction protocol and read in the assay. The same concentrations of Arg<sup>8</sup>-Vasopressin were spiked into the kit Assay Buffer 28, without performing the extraction protocol, for comparison purposes. The efficiency of extraction was calculated as the amount returned off the standard curve divided by the amount returned for the Assay Buffer 28 samples x 100 (after normalizing for endogenous levels).

<b>Assay Buffer 28 Extraction Efficiency</b>			
<b>Vasopressin Spike Level</b>	<b>Spiked Extracted (pg/mL)</b>	<b>Spiked Non-Extracted (pg/mL)</b>	<b>Percent Extraction Recovery</b>
High	214	305	70.2%
Medium	100	140	71.7%
Low	46	61	75.3%

<b>Human Plasma EDTA Extraction Efficiency</b>			
<b>Vasopressin Spike Level</b>	<b>Spiked Extracted (pg/mL)</b>	<b>Percent Extraction Recovery</b>	<b>Plasma vs. AB 28 Extraction Efficiency</b>
High	125	40.90%	58.30%
Medium	93	66.50%	92.70%
Low	36	59.20%	78.60%
<b>Human Serum Extraction Efficiency</b>			
<b>Vasopressin Spike Level</b>	<b>Spiked Extracted (pg/mL)</b>	<b>Percent Extraction Recovery</b>	<b>Plasma vs. AB 28 Extraction Efficiency</b>
High	197	64.40%	91.80%
Medium	93	66.70%	93.00%
Low	28	46.10%	61.20%

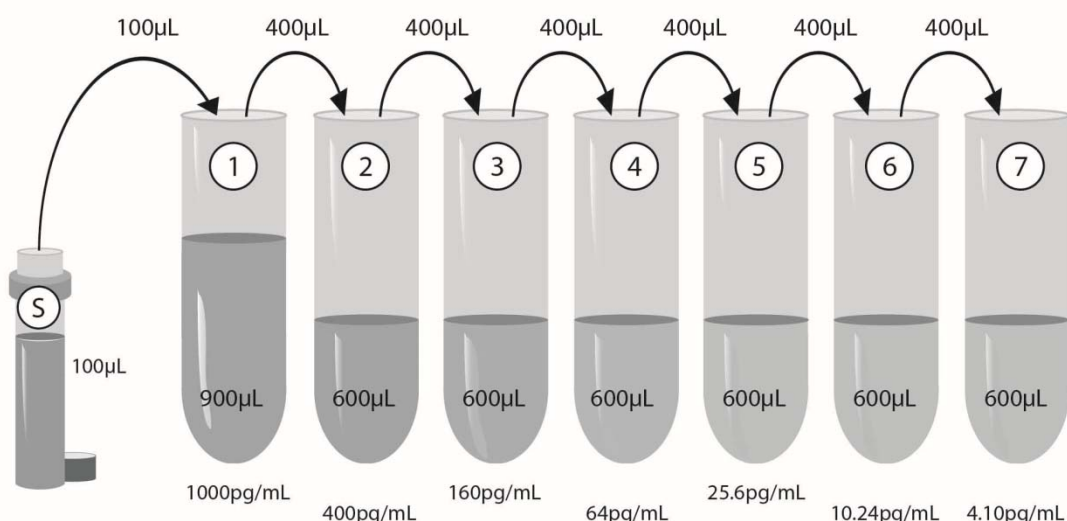
## REAGENT PREPARATION

### 1. Vasopressin Standard

Allow the 10,000 pg/ml Arg<sup>8</sup>-Vasopressin standard solution to warm to room temperature. Label seven 12 x 75mm tubes #1 through #7. Pipet 900µl of Assay Buffer 28 into tube #1 and 600µl into tubes #2 through #7. Add 100µl of the 10,000 pg/ml standard to tube #1. Vortex thoroughly. Add 400µl of tube #1 to tube #2 and vortex thoroughly. Continue this for tubes #3 -#7.



Glass or polypropylene tubes may be used for standard preparation. Avoid polystyrene.



**Diluted standards should be used within 60 minutes of preparation.** The concentrations of Vasopressin in the tubes are labeled above.

### 2. Wash Buffer

Prepare the Wash Buffer by diluting 5ml of the supplied concentrate with 95ml of deionized water. This can be stored at room temperature until the kit expiration date, or for 3 months, whichever is earlier.



Bring all reagents to room temperature for at least 30 minutes prior to opening.



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Pipet standards & samples to the bottom of the wells.



All standards and samples should be run in duplicate.



Prior to the addition of the SA-HRP and substrate, ensure there is no residual wash buffer in the wells. Remaining wash buffer may cause variation in assay results.

## ASSAY PROCEDURE

Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the pouch and seal the ziploc. Store unused wells at 4°C.

1. Add 100µl of Assay Buffer 28 into the NSB and the Bo (0 pg/ml Standard) wells.
2. Add 100µl of Standards #1 through #7 into the appropriate wells.
3. Add 100µl of the Samples into the appropriate wells.
4. Add an additional 50µl of Assay Buffer 28 into the NSB wells.
5. Add 50µl of the blue Conjugate into each well, except the Blank wells.
6. Add 50µl of the yellow Antibody into each well, except the Blank and NSB wells.

**NOTE:** Every well used should be **Green** in color except the NSB wells which should be **Blue**. The Blank wells are empty at this point and have no color.

7. Tap the plate gently to mix. Seal the plate and incubate at 4°C for 18-24 hours.
8. Empty the contents of the plate and wash by adding full well volume (~400 µl) of wash solution to every well. Repeat the wash 2 more times for a total of 3 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
9. Add 200µl of the SA-HRP to every well, except the Blank wells.
10. Seal the plate and incubate at room temperature on a plate shaker for 30 minutes at ~500rpm\*. Wash as above (step 8).
11. Add 200µl of the TMB Substrate to every well.
12. Seal the plate and incubate at room temperature on a plate shaker for 30 minutes at ~500rpm\*.
13. Add 100µl of Stop Solution 2 to every well. This stops the reaction and the plate should be read immediately.
14. Blank the plate reader against the Blank wells, and read the optical density at **450nm**. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all readings.

\* The plate shaker speed was based on a BellCo Mini Orbital Shaker (mod no. 7744-08096). The actual speed of the plate shaker should be such

that the liquid in the plate wells mixes thoroughly, but does not splash out of the well.

## CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of Vasopressin in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic curve fitting program (such as AssayBlaster™, catalog number ADI-28-0002). If data reduction software is not readily available, the concentration of Vasopressin can be calculated as follows:

1. Calculate the average net optical density (OD) bound for each standard and sample by subtracting the average NSB OD from the average OD bound:

$$\text{Average Net OD} = \text{Average Bound OD} - \text{Average NSB OD}$$

2. Plot the Net OD versus the Concentration of Arg<sup>8</sup>-Vasopressin for the standards. Sample concentrations of Arg<sup>8</sup>-Vasopressin may be calculated by interpolation off the standard curve using Net OD values.
3. Alternatively, calculate the binding of each pair of standard wells as a percentage of the maximum binding wells (Bo), using the following formula:

$$\text{Percent Bound} = (\text{Net OD}/\text{Net Bo OD}) \times 100$$

4. Plot Percent Bound versus Concentration of Arg<sup>8</sup>-Vasopressin for the standards. The concentration of Arg<sup>8</sup>-Vasopressin in the unknowns may then also be determined by interpolation off of the binding curve.



Samples with concentrations reading outside of the standard curve range will need to be re-analyzed using a different dilution or more concentrated extract.



Make sure to multiply sample concentrations by the dilution factor used during sample analysis.

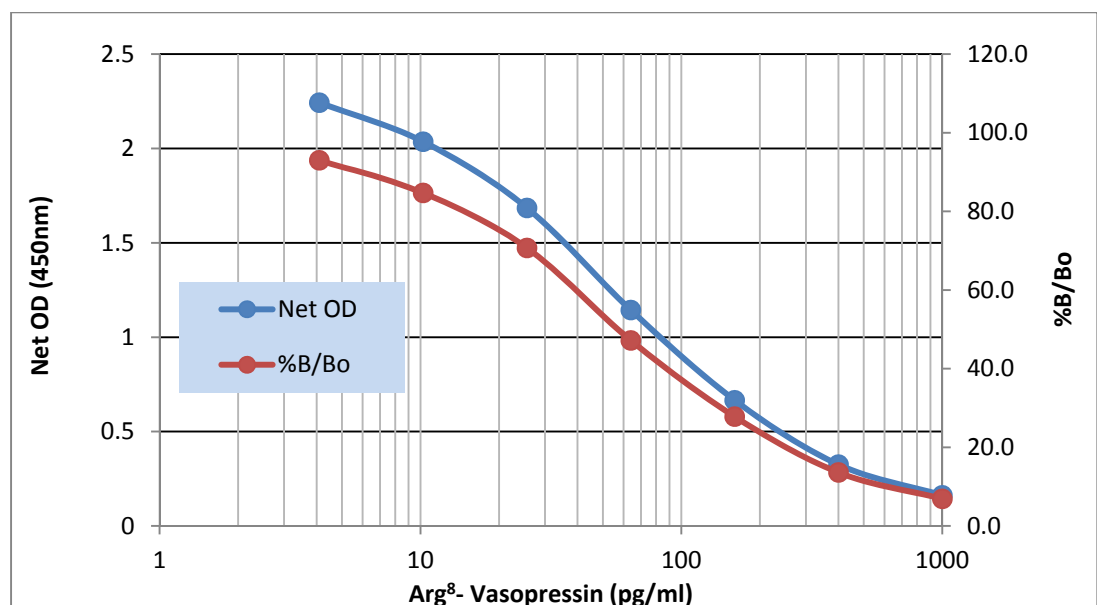
## TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results.

Sample	Arg <sup>8</sup> – Vasopressin (pg/mL)	Net Optical Density (450 nm)	%B/Bo
Bo	0	2.419	---
S1	1000	0.162	6.9
S2	400	0.325	13.6
S3	160	0.666	27.8
S4	64	1.144	47.2
S5	25.6	1.685	70.7
S6	10.24	2.032	84.7
S7	4.096	2.243	93.0
NSB	---	0.031	---

## TYPICAL STANDARD CURVES

Typical standard curves are shown below. These curves must not be used to calculate Arg<sup>8</sup> - Vasopressin concentrations; each user must run a standard curve for each assay.



## PERFORMANCE CHARACTERISTICS

### Sensitivity

The sensitivity of the assay was determined by interpolation from the average of 9 separate standard curves run with replicate data points at each concentration. The sensitivity was determined at 2 standard deviations below the average net OD of 54 zero standard replicates (6 per standard curve). The sensitivity (limit of detection) of the assay is 2.84 pg/ml.

### Linearity

A buffer sample containing Vasopressin was serially diluted 5 times 1:2 in the kit Assay Buffer and measured in the assay. The results follow:

Dilution	Expected (pg/ml)	Observed (pg/ml)	Recovery (%)
Neat	---	390.8	---
1:2	195.4	194.6	100%
1:4	97.7	105.1	108%
1:8	48.8	67.0	137%
1:16	24.4	27.9	114%
1:32	12.2	12.1	99%

The data was plotted graphically as expected Vasopressin concentration versus observed Vasopressin concentration. The line obtained had a slope of 1.0162 with a correlation coefficient of 0.9895.

### Precision

Intra-assay precision was determined by taking samples containing low, medium and high concentrations of Vasopressin and running these samples multiple times (n=22) in the same assay.

The precision numbers listed below represent the percent coefficient of variation for the concentrations of Vasopressin determined in these assays as calculated by a 4 parameter logistic curve fitting program.

Intra-assay precision	
pg/mL	%CV
143.7	6.0%
70.7	6.7%
32.1	14.3%

Inter-assay precision was determined by measuring three samples with low, medium and high concentrations in multiple assays (n=13) over several days.

Inter-assay precision	
pg/mL	%CV
136.2	8.6%
66.4	6.4%
33.0	9.5%

### Cross Reactivities

The % cross reactivity for each related compound was determined by running serial dilutions of each compound (10,000 pg/mL – 10 pg/mL) in the assay, fitting the resulting dose response curve to 4PL curve and determining the ED50. The ED50 of the standard curve was then divided by the determined ED50 of the cross-reactant and multiplied by 100.

Analyte	Cross Reactivity
Arg <sup>8</sup> -Vasopressin	100%
Lys <sup>8</sup> -Vasopressin	9.8%
Oxytocin	<0.001%
TRH	<0.001%
VIP	<0.001%
Leu-Enkephalin	<0.001%
Met-Enkephalin	<0.001%
Mesotocin	<0.001%
Cyclo-Somatostatin	<0.001%
Vasotocin	4.8%
Desmopressin	3.1%
Ser <sup>4</sup> , Ile <sup>8</sup> -Oxytocin	<0.001%



For detailed cross-reactivity protocol, see our website.



## SAMPLE RECOVERIES

Please refer to pages 6-8 for Sample Handling recommendations and Standard preparation. Sample recovery dilutions should only be used to remove matrix interference in media such as Tissue Culture Media.

Plasma and serum samples will need to be extracted and reconstituted in Assay Buffer 28. Upon reconstitution, no further dilution is necessary.

Vasopressin concentrations were measured in tissue culture media. Vasopressin was spiked into the undiluted media and measured neat or following dilution with Assay Buffer 28. Control spikes into Assay Buffer 28 were also measured. The following results were obtained:

Sample	% Recovery*	Recommended Dilution*
Tissue Culture Media	116.2%	no dilution needed
Tissue Culture Media with 10% bovine serum	100.0%	1:2
Assay Buffer 28	111.4%	no dilution needed

\*See Sample Handling instructions on page 6-7 for details.

**REFERENCES**

1. Chard, T, in "An Intro. to Radioimmunoassay & Related Tech.", (1990), 4th Ed., Elsevier, Amsterdam.
2. Tijssen, P, in "Practice & Theory of Enz. Immunoassays", (1985), Elsevier, Amsterdam.
3. Cowley, A.W., et. al, Hypertension, (1981), 3, 93.
4. Clark, G., Wood, P., Merrick, L.,and Lincoln, D.W., Nature, (1979), 282, 746.
5. Malvin, R.L., Fed. Proc., (1971), 30, 1383.
6. Lester, M.C., and Nelson, P.B., Neurosurg., (1981), 8, 735.
7. Schrier, R.W. and Goldberg, J.P., Yale J. Biol. Med., (1980), 53, 525.
8. Chabardes, D, et. al., J. Clin. Invest., (1980), 65, 439.
9. Grill, V and Cerasi, E, J. Biol. Chem., (1974), 249, 41961.
10. Haynes, RC, J. Biol. Chem., (1958), 233, 1220.
11. Szentivanyi, A, J. Allergy, (1968), 42, 203.
12. Hamet, P, et. al, Adv. Cycl. Nucl. Res., (1983), 15, 11.
13. Plaut, M, et. al, Adv. Cycl. Nucl. Res., (1983), 12, 161.
14. Exton, JH, Adv. Cycl. Nucl. Res., (1983), 12, 319.
15. National Committee for Clinical Laboratory Standards Evaluation Protocols, SC1, 1989, NCCLS, Villanova, PA, 19085.



# Product Manual

## NOTES

---



# Product Manual

## **GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
info-usa@enzolifesciences.com

## **EUROPE/ASIA**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)