

### Arg8-Vasopressin ELISA Kit ab205928

[3 References](#) [1 Image](#)

#### Overview

**Product name** Arg8-Vasopressin ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
32.1 pg/mL	22			14.3%
70.7 pg/mL	22			6.7%
143.7 pg/mL	22			6%

Inter-assay

Sample	n	Mean	SD	CV%
33.0	13			9.5%
66.4	13			6.4%
136.2	13			8.6%

**Sample type** Cell culture supernatant, Serum, Plasma

**Assay type** Competitive

**Sensitivity** 2.84 pg/ml

**Range** 4.1 pg/ml - 1000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Tissue Culture Media	116.2	% - %

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Species independent

**Product overview** Abcam's Arg<sup>8</sup>-Vasopressin *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay)

kit is designed for the accurate quantitative measurement of Arg<sup>8</sup>-Vasopressin in tissue culture media.

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with a Biotin conjugated- Arg<sup>8</sup>-Vasopressin antigen and a polyclonal rabbit antibody specific to Arg<sup>8</sup>-Vasopressin. After incubation the excess reagents are washed away. pNpp substrate is added and after a short incubation the enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of Arg<sup>8</sup>-Vasopressin captured in the plate.

## Notes

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). 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, and its release into peripheral blood causes effects upon a number of factors, including emotional stress, posture, blood volume, and temperature. 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.

## Platform

Microplate

## Properties

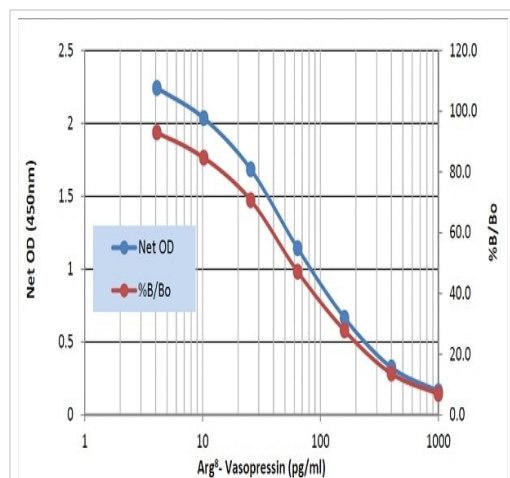
### Storage instructions

Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Assay Buffer	1 x 27ml
Goat anti-rabbit IgG Microplate (12 x 8 wells)	1 unit
Plate Sealer	1 unit
SA-HRP	1 x 20ml
Stop Solution	1 x 10ml
TMB Substrate	1 x 20ml
Vasopressin Biotin Conjugate	1 x 5ml
Vasopressin polyclonal rabbit Antibody	1 x 5ml
Vasopressin Standard	1 x 500µl

## Images

---



Arg8 - Vasopressin Typical Standard Curve

Arg8 - Vasopressin Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors