

# Arg8-Vasopressin

## ELISA kit

Highly sensitive ELISA kit for the detection of Arg<sup>8</sup>-Vasopressin in serum, plasma and tissue culture media.

The Arg<sup>8</sup>-Vasopressin ELISA kit is a colorimetric competitive enzyme immunoassay kit with results overnight + 1 hour. Absorbance is read at 450 nm.

Arginine vasopressin (AVP) is a nonapeptide hormone produced in magnocellular neurons of the hypothalamus, which functions as an antidiuretic in the kidney and systemically as a regulator of vasoconstriction, blood volume, and blood pressure. AVP acts by binding to a set of specific seven transmembrane domain-containing G-protein coupled receptors (GPCR) classified into three groups: V1a (V1), V1b (V3), and V2. V1b receptors stimulate adrenocorticotrophic hormone (ACTH) release in the anterior pituitary, but may also participate in corticotropin-releasing hormone (CRH) release by the hypothalamus. The V1b receptor signals through coupling of Gq/11 alpha subunits, activating phospholipase C and release of intracellular Ca<sup>2+</sup> via phosphoinositide second messengers. V1b receptor overexpression is a common marker for ACTH-secreting corticotrophic adenomas. V2 receptors in the collecting ducts of the kidneys stimulate adenylyl cyclase activity, regulating the reabsorption of water via aquaporin channels.

Citations: 88

[View Online »](#)

### Ordering Information

[Order Online »](#)

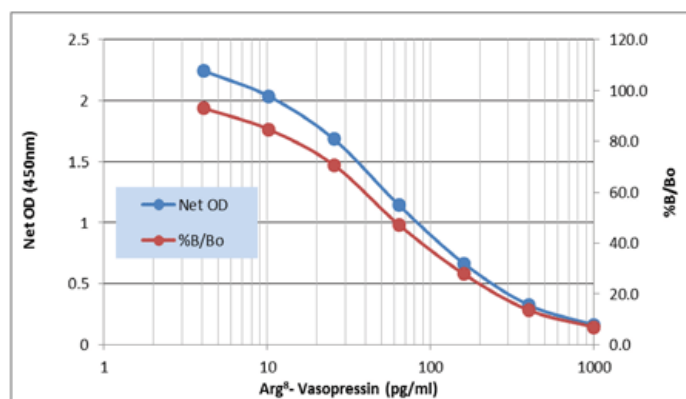
ADI-900-017A

96 wells

### Manuals, SDS & CofA

[View Online »](#)

- High sensitivity, detecting as little as 2.84 pg/ml
- Highly specific with negligible reactivity with oxytocin
- Simple and easy-to-use liquid color-coded reagents reduces error
- Fully quantitative results that surpass semi-quantitative Western blot analysis



## Handling & Storage

**Use/Stability** Store all components at +4°C, except standard at -20°C.

**Shipping** Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Alternative Name** AVP

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of Arg<sup>8</sup>-Vasopressin in serum, plasma, and tissue culture media.

**Assay Time** Overnight + 1 hour

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** GxR IgG Microtiter plate, Conjugate, Antibody, Assay buffer, Wash buffer concentrate, Standard, SA-HRP, TMB Substrate, Stop solution

**Sensitivity** 2.84 pg/ml (range 4.10 – 1,000 pg/ml)

**Species Reactivity** Species independent

**Technical Info / Product Notes**

**Cited samples:**  
[Hormone ELISAs: Cited Samples](#)  
[Wildlife Endocrinology: Cited Sample Types](#)

**UniProt ID** P01185 (human)

**Wavelength** 450 nm

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)