

# Vasopressin receptor

## V<sub>2</sub> polyclonal antibody

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) receptors regulate vasoconstriction, and are expressed in vascular smooth muscle, kidney, myometrium, bladder, adipocytes, hepatocytes, platelets and testis; V1b (V3) receptors stimulate adrenocorticotrophic hormone (ACTH) release in the anterior pituitary, but may also participate in corticotropin-releasing hormone (CRH) release by the hypothalamus; and V2 receptors of the kidney regulate the antidiuretic effects of AVP. 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.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

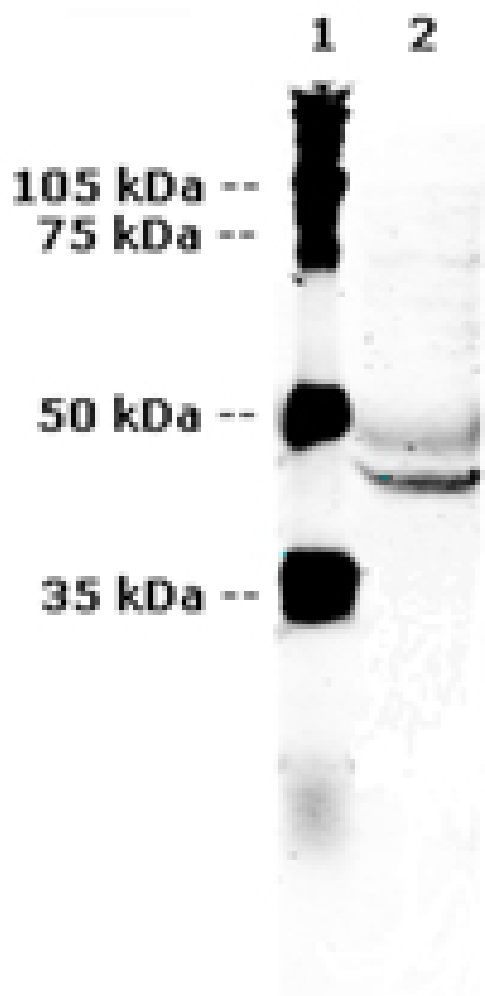
Ordering Information

[Order Online »](#)

ADI-905-812-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW Marker (1) and 20 $\mu$ g of rat kidney extract (2), probed with Vasopressin Receptor (V2) pAb at 2 $\mu$ g/ml.

## Handling & Storage

**Handling** Keep on ice at all times.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** AVPR2

**Application** ELISA, IHC, WB

**Application Notes** Suitable for cell-based and membrane-based ELISA. Predicted MW of ~41kDa. Higher MW species (~50kDa) also observed by Western blot which reflects post-translational modification of the receptor.

**Formulation** Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the N-terminus of rat vasopressin receptor V2.

**Purity Detail** Peptide affinity purified.

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** Q00788

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

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-  
eu@enzolifesciences.com](mailto:info-eu@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)