

Vasopressin receptor

V_{1a} 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 Ca²⁺ 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 »](#)

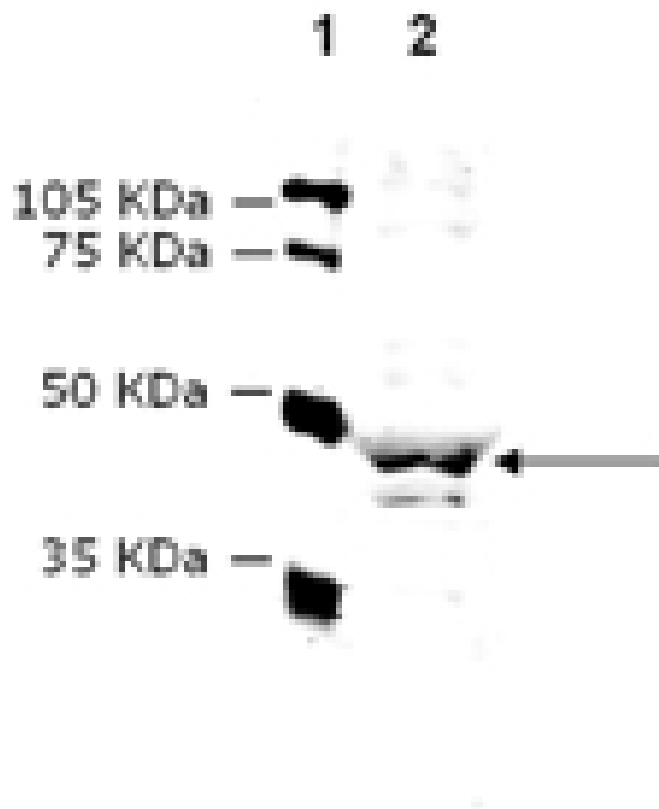
Ordering Information

[Order Online »](#)

ADI-905-811-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: MW Marker (1) and 20µg of rat kidney extract (2), probed with Vasopressin Receptor (V1a) pAb at 1µg/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 AVPR1A

Application WB

Application Notes Predicted MW of ~48kDa. Higher MW species (~50kDa) also observed by Western blot which reflects post-translational modification of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence homology: rhesus macaque, dog, bovine.

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 V1a.

Purity Detail Peptide affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID P30560

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



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)