Arg8-Vasopressin ELISA kit

Highly sensitive ELISA kit for the detection of Arg⁸-Vasopressin in serum, plasma and tissue culture media.

The Arg⁸-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 adrenocorticotropic 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 Ca²⁺ via phosphoinositide second messengers. V1b receptor overexpression is a common marker for ACTH-secreting corticotropic adenomas. V2 receptors in the collecting ducts of the kidneys stimulate adenylyl cyclase activity, regulating the reabsorption of water via aquaporin channels.

Citations: 87

View Online »

Ordering Information

Order Online »

ADI-901-017A

5x96 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 Arg⁸-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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com