

# Angiotensin II ELISA

## kit

Highly sensitive ELISA, detecting as little as 4.6 pg/ml

The Angiotensin II EIA kit is a colorimetric, competitive immunoassay kit with results in 3.5 hours.

Angiotensins are small peptides derived from angiotensinogen. Several of the known angiotensins are established endocrine effectors in the regulation of blood pressure, but they are also known to have other functions locally, in several organs and tissues (paracrine) and at the cellular level (autocrine / intracrine). Angiotensin I (DRVYIHPFHL) has no known effector function but it is an immediate precursor of Angiotensin II (DRVYIHPF). Angiotensin II binds AT1 receptors, which promote vasoconstriction, sodium retention, release of aldosterone, release of Arg-vasopressin, cell proliferation, inflammation, fibrosis, anxiety, and cardiac hypertrophy. Angiotensin A (ARVYIHPF), and Angiotensin III (RVYIHPF), also bind AT1 receptors. Angiotensin (1-7) (DRVYIHP) binds a different receptor called MAS-1 which has opposite effects (vasodilation, natriuresis, antiproliferation, NO release, PGE release, and apoptosis). Angiotensin IV (VYIHPF) binds yet another receptor called AT4 (IRAP), which promotes increase of blood flow, angiogenesis, and natriuresis, and which has also been implicated in memory formation and in the pathogenesis of Alzheimer's disease. The peptide LVV-hemorphin-7 (LVVYPWTQRF), which is not an angiotensin, also binds the AT4 receptor. There are other angiotensin peptides that have been identified, including Ang (1-9) (DRVYIHPFH), Ang (1-12) (DRVYIHPFHLVI), Ang V(3-7) (VYIHP), as well as several other shorter peptides that have undetermined functions.

Citations: 65

[View Online »](#)

## Ordering Information

[Order Online »](#)

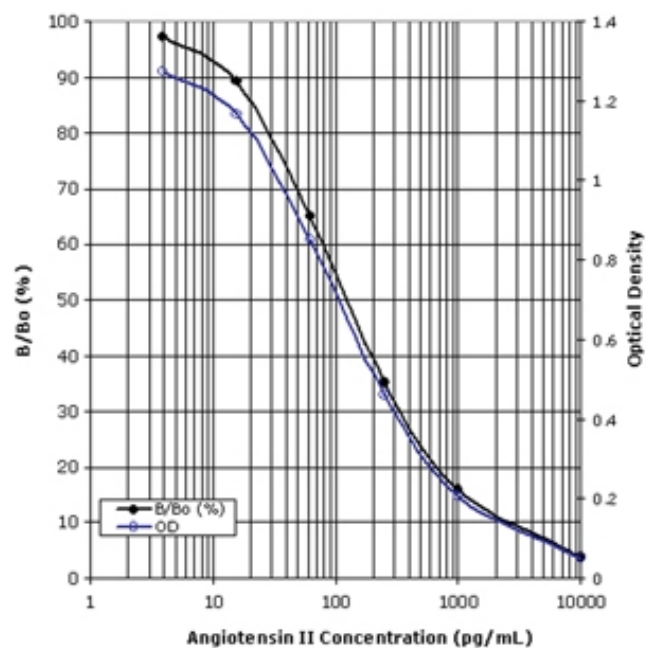
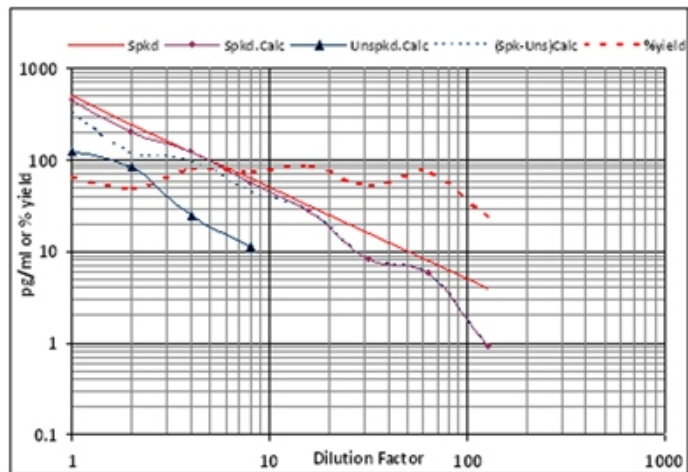
ADI-900-204

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Highly specific with low cross-reactivity to similar peptides
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Easy-to-use liquid color-coded reagents reduce errors
- Ready-to-use reagents and pre-coated plates reduce errors and save your time



## 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

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of species independent Angiotensin II in plasma and serum samples.

**Assay Time** 3.5 hours

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

**Contents** Microtiter Plate, Conjugate, Antibody, Assay Buffer 16, Wash Buffer Concentrate, Standard, TMB Substrate, Stop Solution 2, Streptavidin-HRP

**Crossreactivity** Angiotensin II (100%), Angiotensin A (100%), Angiotensin III (100%), Angiotensin IV (100%), Angiotensin I (0.319%), Angiotensin (1-9) (0.103%), LVV-hemorphin 7 (0.067%), Angiotensin (1-7) (0.053%), Bradykinin (0.022%)

**Quantity** 1 x 96-well plate

**Sensitivity** 4.6 pg/ml (range 3.9-10,000 pg/ml)

**Species Reactivity** Species independent

**Technical Info / Product Notes** **Cited samples:**  
[Wildlife Endocrinology: Cited Sample Types](#)

**UniProt ID** P01019 (human)

### Wavelength



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

### 450 nm

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)

**Last modified: May 29, 2024**

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.