

Angiotensin I ELISA kit

The Angiotensin I ELISA kit is a colorimetric, competitive immunoassay kit

The Angiotensin I 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: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

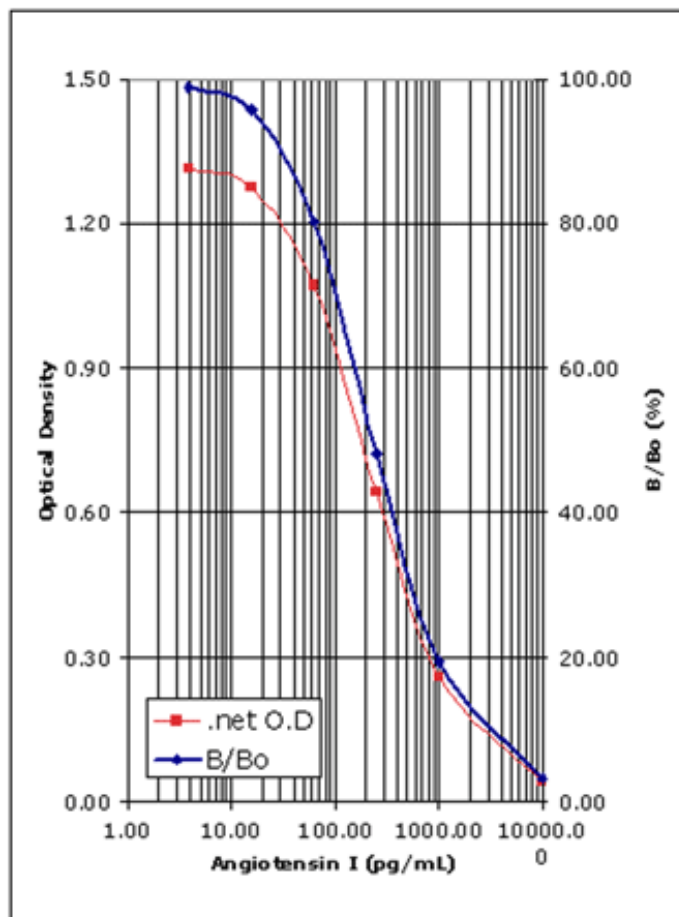
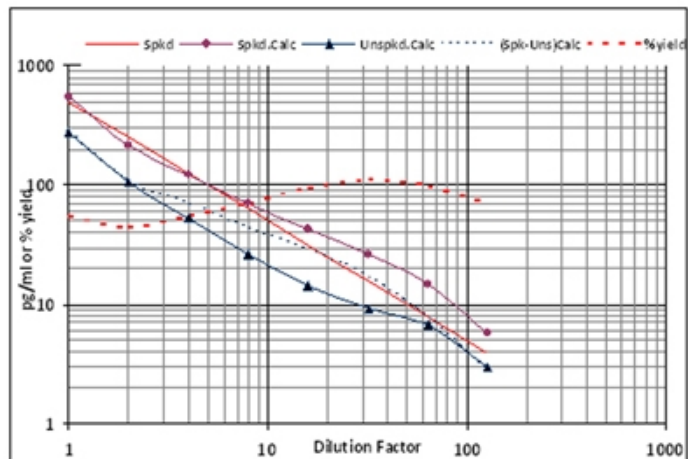
ADI-900-203

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Specific - low cross-reactivity to similar peptides
- Non-radioactive - enzymatic assay format is safe and reliable
- 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	All components of this kit, except the Standard (80□2361) and Conjugate (80□2363), should be stored at +4°C upon receipt.
Short Term Storage	-20°C
Long Term Storage	-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 I 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 I (100%), Angiotensin A (0.182%), Angiotensin II (0.083%), Angiotensin (1-12) (0.025%). Unaffected by Angiotensin (1-9), Angiotensin (1-7), Angiotensin III, Angiotensin IV, LVV-hemorphin 7, Bradykinin.
Quantity	1 x 96-well plate
Sensitivity	4.3 pg/ml (range 3.9-10,000 pg/ml)
Species Reactivity	Species independent
UniProt ID	P01019 (human)
Wavelength	450 nm



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)

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)