

Product datasheet

Angiotensin A ELISA Kit ab136935

[2 Images](#)

Overview

Product name Angiotensin A ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Buffer	20	1297.9pg/ml		7.5%
Buffer	20	655.4pg/ml		4.3%
Buffer	20	55.2pg/ml		10.3%

Inter-assay

Sample	n	Mean	SD	CV%
Buffer		1248.2pg/ml		10.9%
Buffer		612.2pg/ml		5.2%
Buffer		55.6pg/ml		17.7%

Sample type Serum, Plasma

Assay type Competitive

Sensitivity 8.3 pg/ml

Range 9.8 pg/ml - 10000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	94	50pg/ml - 5000pg/ml
Plasma	109	50pg/ml - 5000pg/ml

Assay time 3h 00m

Assay duration

Multiple steps standard assay

Species reactivity**Reacts with:** Human**Product overview**

Abcam's Angiotensin A *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Angiotensin A in plasma and serum.

A goat anti-rabbit IgG polyclonal antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with a solution of Angiotensin A conjugated to biotin, followed by a solution of polyclonal antibody to Angiotensin A. The plate is washed to remove unbound reagents. A solution of streptavidin-HRP conjugate is then added. After further incubation the excess reagents are washed away and TMB substrate is added, which is catalyzed by HRP to generate a yellow color. A stop solution changes this color from yellow to blue, and the intensity of this blue coloration is inversely proportional to the amount of Angiotensin A captured in the plate.

Platform

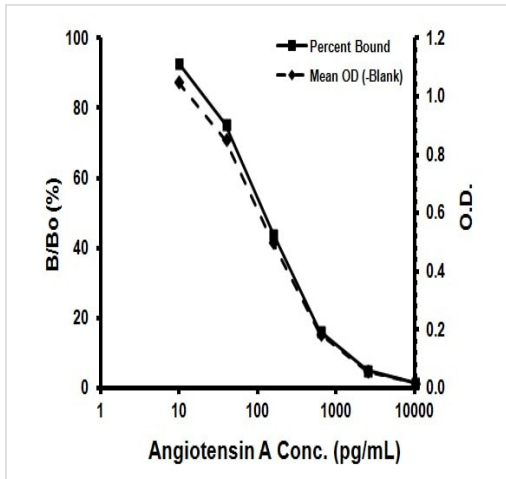
Microplate

Properties**Storage instructions**

Please refer to protocols.

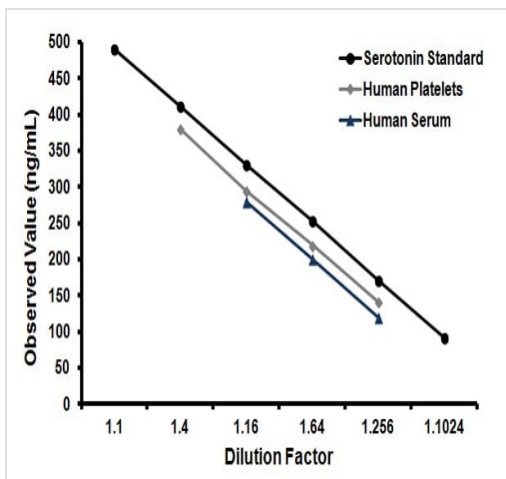
Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Angiotensin A Antibody	1 vial
Angiotensin A Biotin Conjugate	1 vial
Angiotensin A Standard	2 x 40ng
Assay Buffer 16	1 x 30ml
Goat anti-rabbit IgG Microplate (12 x 8 wells)	1 unit
HRP- Streptavidin Conjugate	1 x 12.5µg
Plate Sealer	2 units
Stop Solution 2	1 x 10ml
TMB Substrate	2 x 10ml

Images



Representative standard curve using ab136935.

Typical Standard Curve



Dose-response curves from human plasma and serum diluted into assay buffer were compared to the Angiotensin A standard curve. The parallel response indicates the standard effectively mimics the native protein.

Typical Parallelism Curve

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors