

Product datasheet

Angiotensin I ELISA kit ab136934

1 Image

Overview

Product name Angiotensin I ELISA kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
AGT	20	163.6pg/ml		8.3%
AGT	20	53.2pg/ml		7.3%
AGT	20	25.1pg/ml		10%

Inter-assay

Sample	n	Mean	SD	CV%
AGT		182.2pg/ml		7.9%
AGT		48.1pg/ml		10%
AGT		24.4pg/ml		7.6%

Sample type Serum, Plasma

Assay type Competitive

Sensitivity 4.3 pg/ml

Range 3.9 pg/ml - 10000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Spike	77.3	38pg/ml - 3286pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

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

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with a solution of LVV Hemorphin 7 conjugated to biotin, followed by a solution of polyclonal antibody to LVV Hemorphin 7. 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 LVV Hemorphin 7 captured in the plate.

Notes

Analyte	Sequence	Percent cross reactivities in the range of 0.1 pM - 500 nM
Ang(1-12)	DRVYIHPFHLVI	0.025
Ang I	DRVYIHPFHL	100
Ang(1-9)	DRVYIHPFH	0.000
Ang II	DRVYIHPF	0.083
Ang(1-7)	DRVYIHP	0.000
Ang A	ARVYIHPF	0.182
Ang III	RVYIHPF	0.000
Ang IV	VYIHPF	0.000
LVV-hemorphin 7	LVVYPWTQRF	0.000
Bradykinin	RPPGFSPFR	0.000

Platform

Microplate

Properties

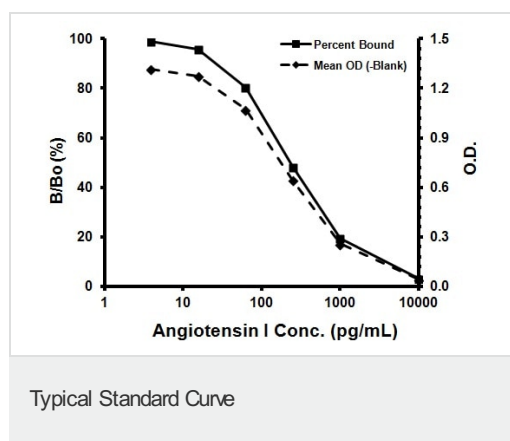
Storage instructions

Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Angiotensin I Biotinylated Conjugate	1 vial
Angiotensin I Polyclonal Antibody	1 vial

Components	1 x 96 tests
Angiotensin I Standard	1 x 1µg
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	1 x 10ml

Images



Representative standard curve using ab136934.

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