

Aldosterone ELISA kit ab136933

★★★★★ [2 Abreviews](#) [28 References](#) [2 图像](#)

概述

产品名称 Aldosterone ELISA试剂盒

检测方法 Colorimetric

精确度

批次内

样品	n	Mean	SD	CV%
Aldosterone	20	84.8pg/ml		4.5%
Aldosterone	20	27.7pg/ml		4.4%
Aldosterone	20	13.9pg/ml		6.6%

批次间

样品	n	Mean	SD	CV%
Aldosterone	13	72.6pg/ml		10.8%
Aldosterone	13	24.4pg/ml		18%
Aldosterone	13	10.8pg/ml		16.3%

样品类型 Urine, Serum, Plasma

检测类型 Competitive

灵敏度 = 4.7 pg/ml

范围 3.9 pg/ml - 250 pg/ml

回收率

特定样本回收率

样品类型	平均%	范围
Urine	= 98.7	7.5pg/ml - 120pg/ml
Serum	= 69.8	7.5pg/ml - 120pg/ml
Plasma	= 108.3	7.5pg/ml - 120pg/ml

实验步骤

Multiple steps standard assay

种属反应性

与反应: Mouse, Rat, Human, Mammals

产品概述

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

A donkey anti-sheep IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with an alkaline phosphatase (AP) conjugated-Aldosterone antigen and a polyclonal sheep antibody specific to Aldosterone. After incubation the excess reagents are washed away. pNpp substrate is added and after a short incubation the enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of Aldosterone captured in the plate.

说明

Compound	% Cross Reactivity
11-Deoxycorticosterone	0.30%
Progesterone	0.20%
Corticosterone	0.19%
Cortisol	≤ 0.001%
DHT	≤ 0.001%
Estradiol	≤ 0.001%
Testosterone	≤ 0.001%

平台

Microplate

性能

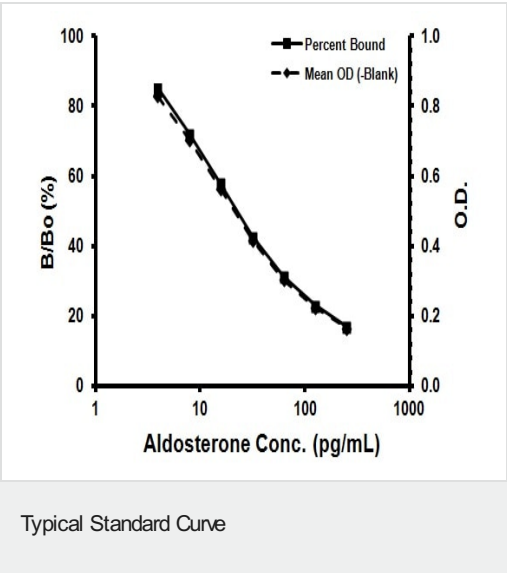
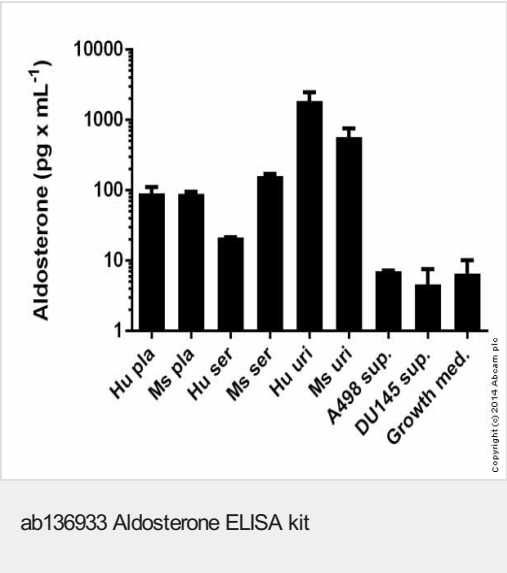
存放说明

Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer Concentrate	1 x 27ml
Aldosterone Alkaline Phosphatase Conjugate	1 x 5ml
Aldosterone Antibody	1 x 5ml
Aldosterone Standard	1 x 250µl
Assay Buffer	1 x 27ml
Donkey anti-sheep IgG Microplate (12 x 8 wells)	1 unit
Plate Sealer	1 unit

组件	1 x 96 tests
pNpp Substrate	1 x 20ml
Stop Solution	1 x 5ml

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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