


Corticosterone ELISA kit ab108821

★★★★★ 3 Abreviews 134 References 2 Images

Overview

Product name	Corticosterone ELISA kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			5.4%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			10.6%
Sample type	Saliva, Milk, Urine, Serum, Plasma			
Assay type	Competitive			
Sensitivity	= 0.28 ng/ml			
Range	0.391 ng/ml - 25 ng/ml			
Recovery	= 98 %			
Assay time	3h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Mouse, Rat, Human, Baboon			
	Predicted to work with: Rabbit, Horse, Cow, Dog, Pig 			
Product overview	Corticosterone ELISA kit is a competitive Enzyme-Linked Immunosorbent Assay designed for the quantitative measurement of Corticosterone levels in plasma, serum, urine, milk, and saliva samples.			

A Corticosterone specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Corticosterone is added and then followed by washing with wash buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is inversely

proportional to the amount of Corticosterone captured in plate.

**Please note the concentrations of the kit components are lot-specific and the end user should always refer to the vial label**

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

#### Platform

Microplate

#### Properties

#### Storage instructions

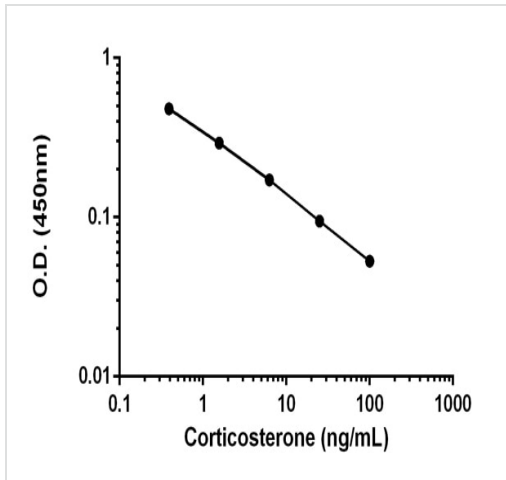
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl	1 x 80µl
10X Diluent S Concentrate	1 x 20ml	1 x 20ml
2X Biotinylated Corticosterone Protein (Lyophilized)	1 vial	1 vial
20X Wash Buffer Concentrate	1 x 30ml	1 x 30ml
Chromogen Substrate	1 x 7ml	1 x 7ml
Corticosterone Microplate (12 x 8 well strips)	1 unit	1 unit
Corticosterone Standard	1 vial	1 vial
Sealing Tapes	3 units	3 units
Stop Solution	1 x 11ml	1 x 11ml

#### Relevance

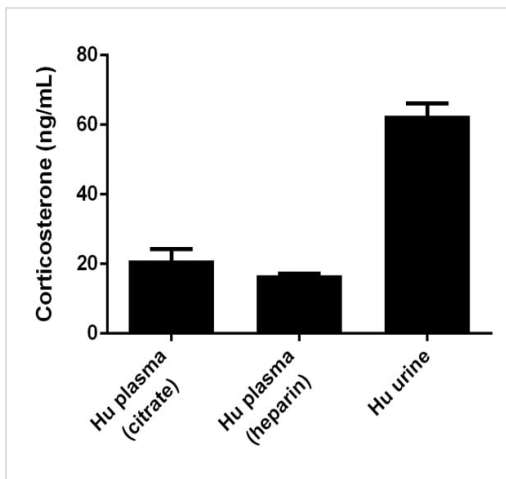
Corticosterone is the adrenal steroid, the major glucocorticoid. Glucocorticoid hormones are also known as corticosteroid hormones and are synthesized mainly in the adrenal cortex; however, more recently the enzymes involved in their synthesis have been found in a variety of cells and tissues, including the heart. The effects of these hormones are mediated via both cytoplasmic mineralocorticoid receptors (MRs) and glucocorticoid receptors (GRs), which act as ligand-inducible transcription factor. Corticosterone has profound effect on the structure and function of the hippocampus. Brain corticosterone action through the glucocorticoid receptor may involve memory storage. Emotional stress might cause increases in plasma corticosterone.

#### Images



Competitive ELISA with ab108821.

Competitive ELISA - Corticosterone (ab108821)



Corticosterone measured in biological fluids showing quantity (ng) per mL of tested sample.

Competitive ELISA - Corticosterone ELISA kit  
(ab108821)

**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