

Corticosterone ELISA kit ab108821

★★★★★ 3 Abreviews 134 References 画像数 2

製品の概要

製品名	Corticosterone ELISA kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				5.4%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				10.6%
サンプルの種類	Saliva, Milk, Urine, Serum, Plasma			
アッセイタイプ	Competitive			
検出感度	= 0.28 ng/ml			
検出範囲	0.391 ng/ml - 25 ng/ml			
添加回収試験	= 98 %			
全工程の試験時間	3h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Mouse, Rat, Human, Baboon			
	交差が予測される動物種: Rabbit, Horse, Cow, Dog, Pig			
製品の概要	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.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

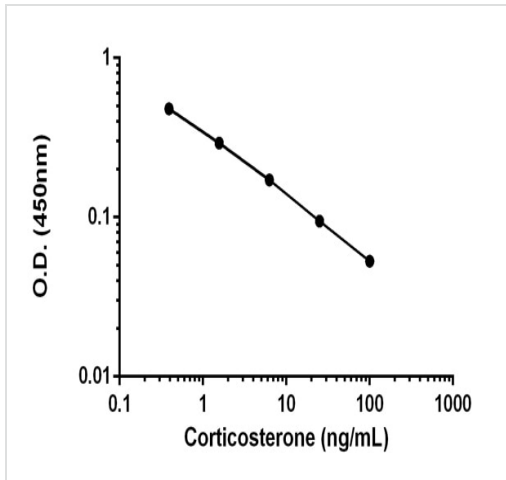
Store at -20°C. Please refer to protocols.

内容	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

#### 関連性

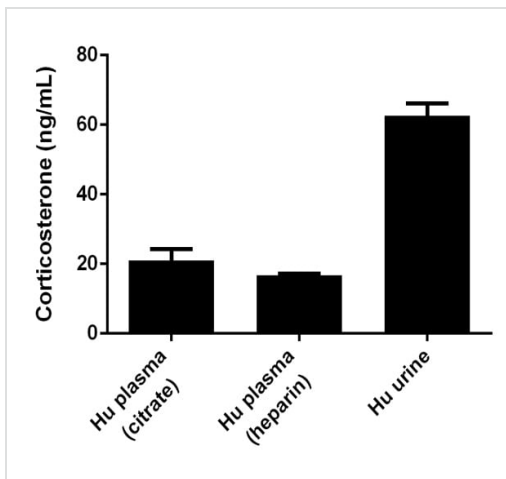
Corticosterone is the adrenal steroid, the major glucocorticoid. Glucocorticoid hormones are also known as corticosteroid hormones and are synthesized mainly in the adrenal cortex; however, more recently the enzymes involved in their synthesis have been found in a variety of cells and tissues, including 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.

#### 画像



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.co.jp/abpromise> or contact our technical team.

#### **Terms and conditions**

---

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