

Product datasheet

Corticosterone ELISA kit ab108821

★ ★ ★ ☆ ☆ 1 Abreviews 23 References 2 Images

Overview

Product name Corticosterone ELISA kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | | | | 5.3% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | 10.6% |

Sample type Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma

Assay type Competitive

Sensitivity = 0.28 ng/ml

Range 2 ng/ml - 22 ng/ml

Recovery = 101 %

Assay time 3h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse, Rat, Human, Baboon

Product overview

Corticosterone *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab108821) is designed for the quantitative measurement of Corticosterone levels in plasma, serum, urine, milk, saliva and cell culture supernatant.

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.

Tested applications**Suitable for:** Competitive ELISA**Platform**

Microplate

Properties**Storage instructions**

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

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

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.

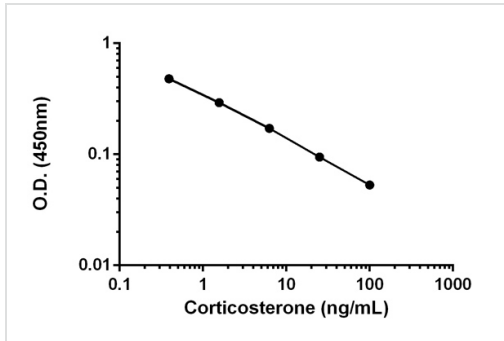
Applications

Our [Abpromise guarantee](#) covers the use of **ab108821** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

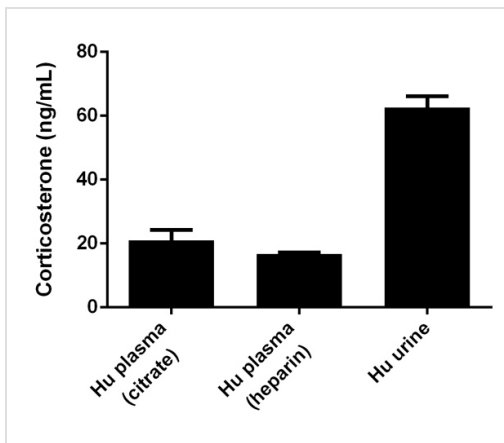
| Application | Abreviews | Notes |
|-------------------|-----------|--|
| Competitive ELISA | | Use at an assay dependent concentration. |

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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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