

Cortisol ELISA Kit ab108674

1 References 1 图像

概述

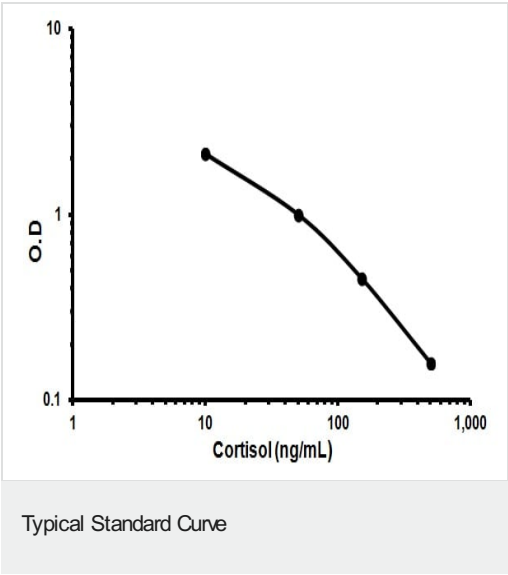
产品名称	Cortisol ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Urine	60		
				CV%
				<=6.5%
	批次间			
	样品	n	Mean	SD
	Urine	30		
				CV%
				<=7.2%
样品类型	Urine			
检测类型	Competitive			
灵敏度	2.95 ng/ml			
范围	10 ng/ml - 500 ng/ml			
回收率	107.48 %			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	<p>Abcam's Cortisol <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Cortisol in urine.</p> <p>A 96-well plate has been precoated with anti-Cortisol. Samples and the Cortisol-HRP conjugate are added to the wells, where Cortisol in the sample competes with the added Cortisol-HRP for antibody binding. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is inversely proportional to the amount of Cortisol in the sample and the intensity is measured at 450 nm.</p>			
平台	Microplate			

性能

存放说明Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Wash Solution	1 x 50ml
Anti-Cortisol IgG Coated Microplate (12 x 8 wells)	1 unit
Cortisol Standard 0 - 0 ng/mL	1 x 4ml
Cortisol Standard 1 - 1 ng/mL	1 x 1ml
Cortisol Standard 2 - 5 ng/mL	1 x 1ml
Cortisol Standard 3 - 30 ng/mL	1 x 1ml
Cortisol Standard 4 – 200 ng/mL	1 x 1ml
Cortisol-HRP Conjugate	1 x 33ml
Cover foils	1 unit
High control	1 x 1ml
Low control	1 x 1ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

图片



Representative Standard Curve using ab108674

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