

Human Complement C4 ELISA Kit ab108824

8 References 2 图像

概述

产品名称	人Complement C4 ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall			5.1%
	批次间			
	样品	n	Mean	SD
	Overall			10.1%
样品类型	Serum, Plasma			
检测类型	Competitive			
灵敏度	= 0.53 µg/ml			
范围	160 µg/ml - 480 µg/ml			
回收率	94 %			
检测时间	3h 00m			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Complement C4 Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab108834) is designed for the quantitative measurement of Complement C4 levels in plasma and serum.			

A Complement C4 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Complement C4 is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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 Complement C4 captured in plate.

Reactivity with C4 cleavage products (C4a, C4b, C4d) has not been determined.

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

平台

Microplate

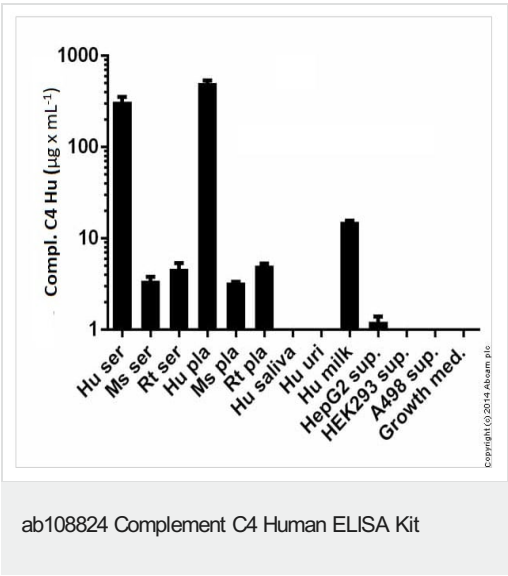
性能

存放说明

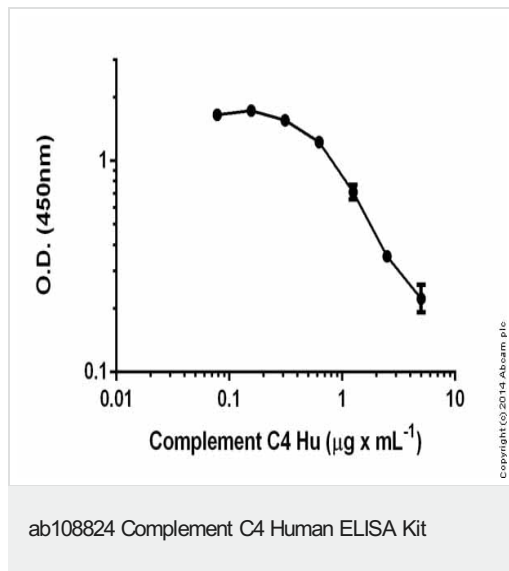
Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	1 x 30ml
2X Biotinylated Human Complement C4 (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
Complement C4 Microplate (12 x 8 well strips)	1 unit
Complement C4 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

图片



Complement C4 measured in biological fluids showing quantity (microgram) per mL of tested sample



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