

★★★★★ 1 Abreviews 5 References 1 图像

概述

产品名称 兔C Reactive蛋白(PTX1) ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 10%

样品类型	Serum, Plasma, Other biological fluids
------	--

检测类型	Sandwich (quantitative)
------	-------------------------

灵敏度 2.0828 ng/ml

范围	6.25 ng/ml - 200 ng/ml
----	------------------------

回收率	特定样本回收率
-----	---------

样品类型	平均%	范围
Serum	> 85	% - %

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rabbit

产品概述 Abcam's C Reactive Protein (PTX1) Rabbit ELISA Kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring C Reactive Protein levels in rabbit serum and plasma.

In this assay the C Reactive Protein present in samples reacts with the anti-C Reactive Protein antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-C Reactive Protein antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound C Reactive Protein. Following another washing step, the enzyme bound

to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of C Reactive Protein in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of C Reactive Protein in the test sample. The quantity of C Reactive Protein in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

平台 Microplate

性能

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

组件	1 x 96 tests
100X HRP-conjugated anti-rabbit C Reactive Protein antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Rabbit C Reactive Protein Calibrator (lyophilized)	1 vial
Rabbit C Reactive Protein ELISA Microplate	1 unit
Stop Solution	1 x 12ml

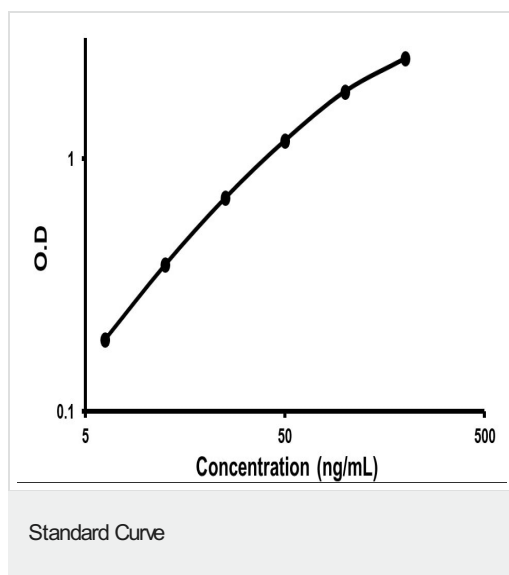
功能 Displays several functions associated with host defense: it promotes agglutination, bacterial capsular swelling, phagocytosis and complement fixation through its calcium-dependent binding to phosphorylcholine. Can interact with DNA and histones and may scavenge nuclear material released from damaged circulating cells.

组织特异性 Found in plasma.

序列相似性 Belongs to the pentaxin family.
Contains 1 pentaxin domain.

细胞定位 Secreted.

图片



Representative standard curve using ab157726 C Reactive Protein Rabbit ELISA Kit

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