

Rabbit C Reactive Protein (PTX1) ELISA Kit ab157726

★★★★★ 1 Abreviews 5 References 画像数 1

医薬用外毒物

医薬用外劇物

製品の概要

製品名	Rabbit C Reactive Protein (PTX1) 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	CV%
	Overall				< 10%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Overall				< 10%
サンプルの種類	Serum, Plasma, Other biological fluids				
アッセイタイプ	Sandwich (quantitative)				
検出感度	2.0828 ng/ml				
検出範囲	6.25 ng/ml - 200 ng/ml				
添加回収試験	特定サンプルでの回収試験				
	サンプルの種類		平均 %	測定範囲	
	Serum		> 85	% - %	
ステップ	Multiple steps standard assay				
種交差性	交差種: Rabbit				
製品の概要	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.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	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

機能

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.

組織特異性

Found in plasma.

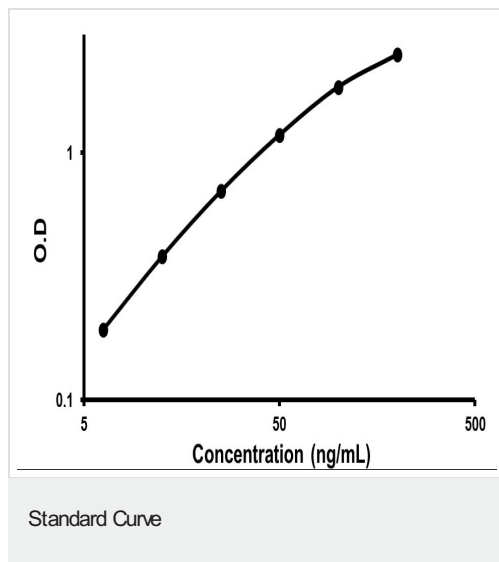
配列類似性

Belongs to the pentaxin family.
Contains 1 pentaxin domain.

細胞内局在

Secreted.

画像



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.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