

Pig CRP ELISA Kit ab205089

2 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	Pig CRP 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%
	Serum				< 10%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Serum				< 10%
サンプルの種類	Serum, Hep Plasma, EDTA Plasma, Cit plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	2.15 ng/ml				
検出範囲	6.25 ng/ml - 200 ng/ml				
添加回収試験	> 85 %				
	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		6.25ng/ml - 200ng/ml	
全工程の試験時間	1h 10m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Pig				
製品の概要	CRP in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CRP protein in pig biological samples.				
特記事項	In this assay the CRP present in samples reacts with the anti-CRP antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-CRP antibodies conjugated with horseradish peroxidase (HRP), are added. These				

enzyme-labeled antibodies form complexes with the previously bound CRP. 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 CRP in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of CRP in the test sample. The quantity of CRP 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

Pre-coated microplate (12 x 8 well strips)

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 Enzyme Antibody	1 x 150µl
1X Diluent Concentration	1 x 60ml
20X Wash Buffer Concentrate	1 x 50ml
Anti-Pig CRP ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Pig CRP Calibrator (Lyophilized)	1 vial
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

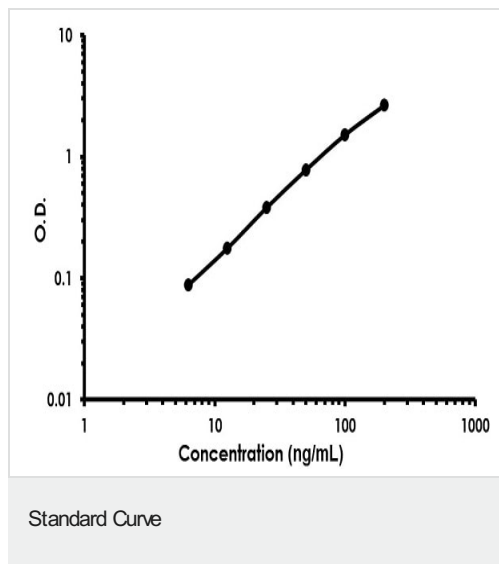
Belongs to the pentaxin family.

Contains 1 pentaxin domain.

細胞内局在

Secreted.

画像



Representative standard curve using ab205089 Pig CRP 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