

★★★★★ **1 Abreviews** **5 References** **1 Image**

Overview

| | |
|---------------------|--|
| Product name | Rabbit C Reactive Protein (PTX1) ELISA Kit |
|---------------------|--|

| | |
|------------------|--------------|
| Detection method | Colorimetric |
|------------------|--------------|

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

| | |
|--------------------|--|
| Sample type | Serum, Plasma, Other biological fluids |
|--------------------|--|

| | |
|-------------------|-------------------------|
| Assay type | Sandwich (quantitative) |
|-------------------|-------------------------|

Sensitivity 2.0828 ng/ml

| | |
|--------------|------------------------|
| Range | 6.25 ng/ml - 200 ng/ml |
|--------------|------------------------|

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|-------------|-----------|-------|
| Serum | > 85 | % - % |

| | |
|-----------------------|-------------------------------|
| Assay duration | Multiple steps standard assay |
|-----------------------|-------------------------------|

Species reactivity **Reacts with:** Rabbit

Product overview 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.

1

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.

Platform Microplate

Properties

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

| Components | 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 |

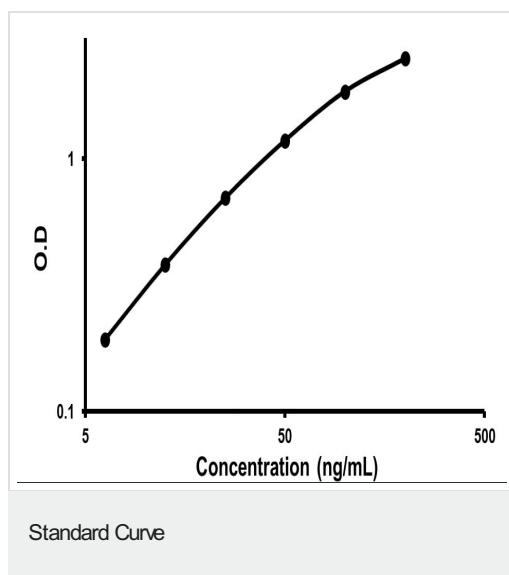
Function Displays several functions associated with host defense: it promotes agglutination, bacterial capsular swelling, phagocytosis and complement fixation through its calcium-dependent binding to phosphorylcholine. Can interact with DNA and histones and may scavenge nuclear material released from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

Tissue specificity Found in plasma.

Sequence similarities Belongs to the pentaxin family.
Contains 1 pentaxin domain.

Cellular localization Secreted.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors