

# Mouse C Reactive Protein ELISA Kit (PTX1) ab157712

[12 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Mouse C Reactive Protein ELISA Kit (PTX1)			
<b>Detection method</b>	Colorimetric			
<b>Precision</b>	Intra-assay			
	<b>Sample</b>	<b>n</b>	<b>Mean</b>	<b>SD</b>
	Overall			< 10%
	Inter-assay			
	<b>Sample</b>	<b>n</b>	<b>Mean</b>	<b>SD</b>
	Overall			< 10%
<b>Sample type</b>	Urine, Serum, Plasma			
<b>Assay type</b>	Sandwich (quantitative)			
<b>Sensitivity</b>	0.1691 ng/ml			
<b>Range</b>	0.78 ng/ml - 25 ng/ml			
<b>Recovery</b>	Sample specific recovery			
	<b>Sample type</b>	<b>Average %</b>	<b>Range</b>	
	Serum	> 85	% - %	
<b>Assay duration</b>	Multiple steps standard assay			
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse			
<b>Product overview</b>	Abcam's C Reactive Protein (PTX1) Mouse ELISA kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring C Reactive Protein in mouse biological samples.			

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.

**Platform** Microplate

## Properties

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

Components	1 x 96 tests
100X HRP-conjugated anti-mouse 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
Mouse C Reactive Protein Calibrator (Lyophilized)	1 vial
Mouse C Reactive Protein ELISA Microplate	1 unit
Stop Solution	1 x 12ml

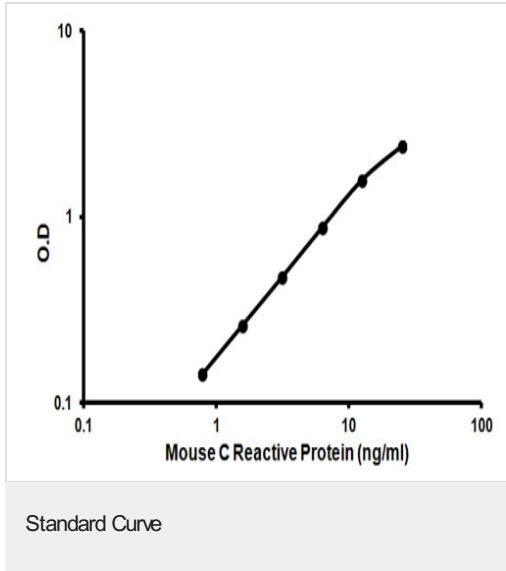
**Function** Displays several functions associated with host defense: it promotes agglutination, bacterial capsular swelling, phagocytosis and complement fixation through its calcium-dependent binding to phosphorylcholine. Can interact with DNA and histones and may scavenge nuclear material released 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 ab157712 C Reactive Protein  
Mouse 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