

### Rat C Reactive Protein ELISA Kit (PTX1) ab108827

[18 References](#) [画像数 1](#)

#### 製品の概要

**製品名** Rat C Reactive Protein ELISA Kit (PTX1)

**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 |   |     |    | 3.4% |

Inter-Assay (日差再現性)

| サンプル    | N | 平均値 | SD | CV%  |
|---------|---|-----|----|------|
| Overall |   |     |    | 8.5% |

**サンプルの種類** Cell culture supernatant, Urine, Serum, Plasma

**アッセイタイプ** Sandwich (quantitative)

**検出感度** = 0.25 ng/ml

**検出範囲** 0.391 ng/ml - 25 ng/ml

**添加回収試験** 95 %

**全工程の試験時間** 3h 30m

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Rat

**製品の概要** Abcam's C Reactive Protein (PTX1) Rat *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of C Reactive Protein in urine, plasma, serum, and cell culture supernatants.

A C Reactive Protein specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a C Reactive Protein specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of C Reactive Protein captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

試験プラットフォーム

Microplate

## 製品の特性

### 保存方法

Store at -20°C. Please refer to protocols.

| 内容   | 1 x 96 tests |
|--|--------------|
| 100X Streptavidin-Peroxidase Conjugate             | 1 x 80µl     |
| 10X Diluent M Concentrate                          | 1 x 30ml     |
| 20X Wash Buffer Concentrate                        | 2 x 30ml     |
| 50X Biotinylated Rat C Reactive Protein Antibody   | 1 x 120µl    |
| C Reactive Protein Microplate (12 x 8 well strips) | 1 unit       |
| C Reactive Protein Standard                        | 1 vial       |
| Chromogen Substrate                                | 1 x 7ml      |
| Sealing Tapes                                      | 3 units      |
| Stop Solution                                      | 1 x 11ml     |

### 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.

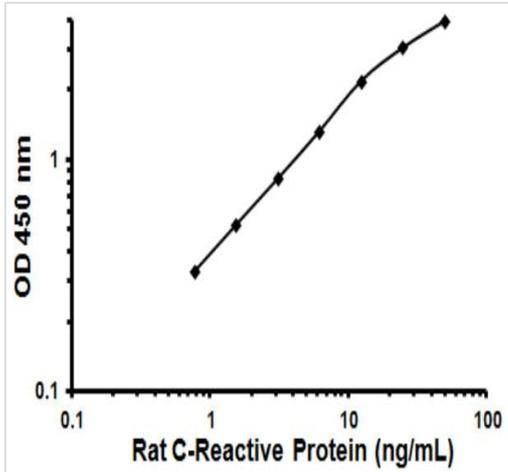
### 配列類似性

Belongs to the pentaxin family.  
Contains 1 pentaxin domain.

### 細胞内局在

Secreted.

## 画像



Representative Standard Curve using ab108827.

Typical Standard Curve

**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