

## Product datasheet

# Human C Reactive Protein ELISA Kit (PTX1) ab108826

1 References [画像数 1](#)

### 製品の概要

<b>製品名</b>	Human C Reactive Protein ELISA Kit (PTX1)																				
<b>検出方法</b>	Colorimetric																				
<b>再現性</b>	<p style="text-align: right;">Intra-Assay (同時再現性)</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>4%</td> </tr> </tbody> </table> <hr/> <p style="text-align: right;">Inter-Assay (日差再現性)</p> <table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>11.3%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				4%	サンプル	N	平均値	SD	CV%	Overall				11.3%
サンプル	N	平均値	SD	CV%																	
Overall				4%																	
サンプル	N	平均値	SD	CV%																	
Overall				11.3%																	
<b>サンプルの種類</b>	Cell culture supernatant, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid																				
<b>アッセイタイプ</b>	Sandwich (quantitative)																				
<b>検出感度</b>	= 0.2 ng/ml																				
<b>検出範囲</b>	0.25 ng/ml - 16 ng/ml																				
<b>添加回収試験</b>	96 %																				
<b>全工程の試験時間</b>	4h 00m																				
<b>ステップ</b>	Multiple steps standard assay																				
<b>種交差性</b>	交差種: Human																				
<b>製品の概要</b>	<p>Abcam's C Reactive Protein (PTX1) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of C Reactive Protein concentrations in Human plasma, serum, saliva, milk, urine and cell culture supernatants.</p> <p>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.</p>																				

Get results in 90 minutes with Human C Reactive Protein ELISA Kit ([ab181416](#)) from our SimpleStep ELISA® range.

#### アプリケーション

**適用あり:** Sandwich ELISA

#### 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 N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human 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 8ml
Sealing Tapes	3 units
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

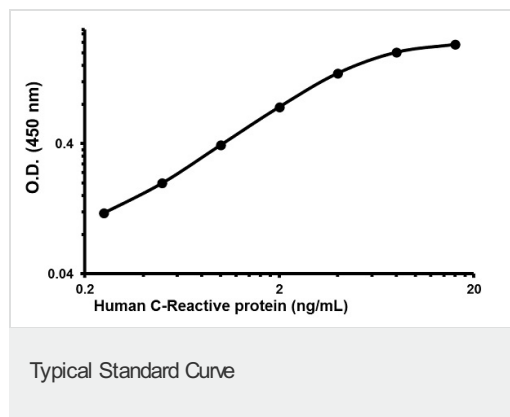
Secreted.

#### アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108826** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.



Representative Standard Curve using ab108826

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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