

Product datasheet

Human C Reactive Protein ELISA Kit (PTX1) ab108826

1 References [画像数 1](#)

製品の概要

製品名	Human C Reactive Protein ELISA Kit (PTX1)																				
検出方法	Colorimetric																				
再現性	<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 style="width: 20%;">サンプル</th> <th style="width: 5%;">N</th> <th style="width: 20%;">平均値</th> <th style="width: 10%;">SD</th> <th style="width: 45%;">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 style="width: 20%;">サンプル</th> <th style="width: 5%;">N</th> <th style="width: 20%;">平均値</th> <th style="width: 10%;">SD</th> <th style="width: 45%;">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%																	
サンプルの種類	Cell culture supernatant, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid																				
アッセイタイプ	Sandwich (quantitative)																				
検出感度	= 0.2 ng/ml																				
検出範囲	0.25 ng/ml - 16 ng/ml																				
添加回収試験	96 %																				
全工程の試験時間	4h 00m																				
ステップ	Multiple steps standard assay																				
種交差性	交差種: Human																				
製品の概要	<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>																				

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.

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

製品の特徴

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

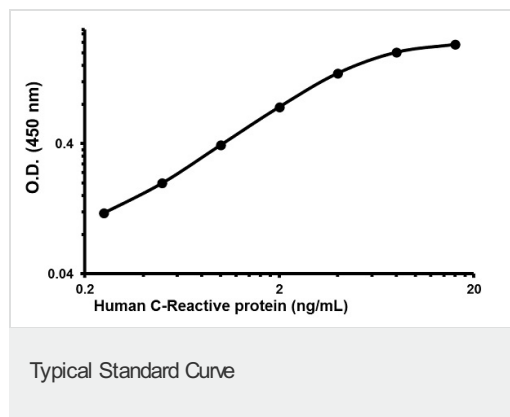
Secreted.

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

Typical Standard Curve

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 <http://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