

### Human Protein C ELISA Kit ab137988

画像数 1

#### 製品の概要

製品名	Human Protein C ELISA Kit										
検出方法	Colorimetric										
再現性	Intra-Assay (同時再現性)										
	<table border="1"> <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>Intra-assay</td> <td></td> <td></td> <td></td> <td>5.9%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Intra-assay				5.9%
サンプル	N	平均値	SD	CV%							
Intra-assay				5.9%							
	Inter-Assay (日差再現性)										
	<table border="1"> <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>Inter-assay</td> <td></td> <td></td> <td></td> <td>10.6%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Inter-assay				10.6%
サンプル	N	平均値	SD	CV%							
Inter-assay				10.6%							
サンプルの種類	Serum, Plasma										
アッセイタイプ	Competitive										
検出感度	55 ng/ml										
検出範囲	0.094 µg/ml - 6 µg/ml										
添加回収試験	97 %										
全工程の試験時間	3h 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 Protein C Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Protein C in plasma and serum.</p> <p>A Protein C specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Protein C is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 inversely proportional to the amount of Protein C captured in plate.</p>										

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
1X Biotinylated Human Protein C (Lyophilized)	1 vial
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 20ml
20X Wash Buffer Concentrate	1 x 30ml
Chromogen Substrate	1 x 7ml
Protein C Microplate (12 x 8 well strips)	1 unit
Protein C Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11 ml

### 機能

Protein C is a vitamin K-dependent serine protease that regulates blood coagulation by inactivating factors Va and VIIIa in the presence of calcium ions and phospholipids (PubMed:25618265). Exerts a protective effect on the endothelial cell barrier function (PubMed:25651845).

### 組織特異性

Plasma; synthesized in the liver.

### 関連疾患

Thrombophilia due to protein C deficiency, autosomal dominant  
Thrombophilia due to protein C deficiency, autosomal recessive

### 配列類似性

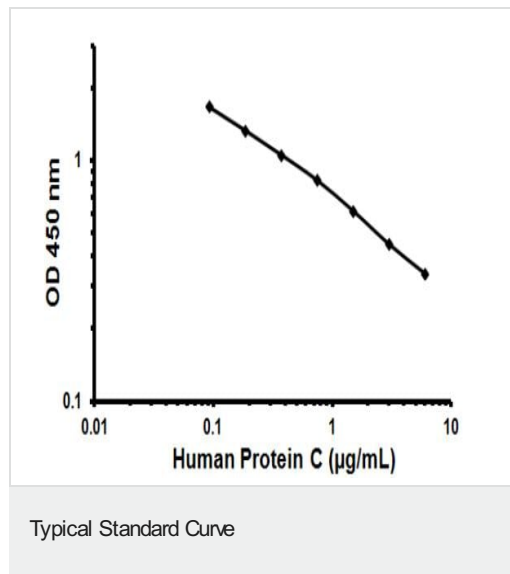
Belongs to the peptidase S1 family.  
Contains 2 EGF-like domains.  
Contains 1 Gla (gamma-carboxy-glutamate) domain.  
Contains 1 peptidase S1 domain.

### 翻訳後修飾

The vitamin K-dependent, enzymatic carboxylation of some Glu residues allows the modified protein to bind calcium.  
N- and O-glycosylated. Partial (70%) N-glycosylation of Asn-371 with an atypical N-X-C site produces a higher molecular weight form referred to as alpha. The lower molecular weight form, not N-glycosylated at Asn-371, is beta. O-glycosylated with core 1 or possibly core 8 glycans. The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.  
May be phosphorylated on a Ser or Thr in a region (AA 25-30) of the propeptide.

### 細胞内局在

Secreted. Golgi apparatus. Endoplasmic reticulum.



Representative Standard Curve using ab137988

**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