

Human Protein C ELISA Kit ab137987

1 References 1 Image

Overview

Product name	Human Protein C ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			= 3.5%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			= 10.2%
Sample type	Cell culture supernatant, Saliva, Milk, Urine, Tissue, Cell Lysate			
Assay type	Sandwich (quantitative)			
Sensitivity	= 2 ng/ml			
Range	12.5 ng/ml - 100 ng/ml			
Recovery	= 97 %			
Assay time	4h 0m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Protein C Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Protein C in urine, saliva, milk, cell culture supernatant, cell lysate and tissue samples.			

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 a Protein C specific biotinylated detection antibody 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 directly proportional to the amount of Protein C captured in plate.

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

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	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 Protein C Antibody	1 x 120µl
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 11ml

**Function** 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).

**Tissue specificity** Plasma; synthesized in the liver.

**Involvement in disease** Thrombophilia due to protein C deficiency, autosomal dominant  
Thrombophilia due to protein C deficiency, autosomal recessive

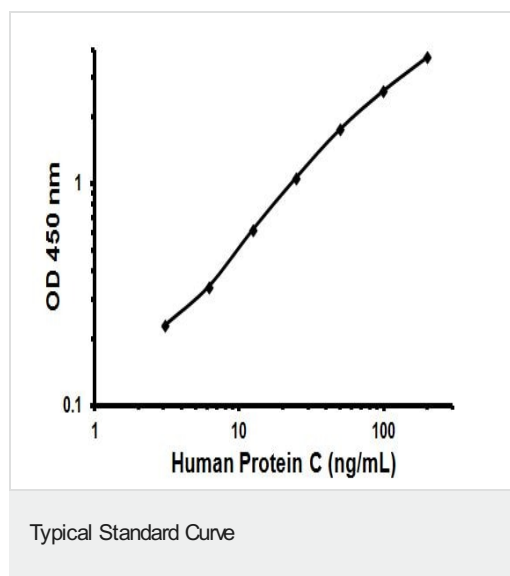
**Sequence similarities** Belongs to the peptidase S1 family.  
Contains 2 EGF-like domains.  
Contains 1 Gla (gamma-carboxy-glutamate) domain.  
Contains 1 peptidase S1 domain.

**Post-translational modifications** 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.

**Cellular localization** Secreted. Golgi apparatus. Endoplasmic reticulum.

## Images

---



Representative Standard Curve using ab137987

**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