

Human Factor X ELISA Kit ab108832

2 References 4 Images

Overview

Product name	Human Factor X ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			3.8%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			8.8%
Sample type	Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	= 1.5 ng/ml			
Range	0.781 ng/ml - 50 ng/ml			
Recovery	97.5 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Human Factor <i>Xin vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor X concentrations in plasma, serum, milk, urine, saliva, and cerebrospinal fluid.			

A Factor X specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor X 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 Factor X 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 Biotinylated Antibody	1 x 60µl
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
Chromogen Substrate	1 x 7ml
Factor X Microplate (12 x 8 well strips)	1 unit
Factor X Standard	1 x 65ng
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Function** Factor Xa is a vitamin K-dependent glycoprotein that converts prothrombin to thrombin in the presence of factor Va, calcium and phospholipid during blood clotting.

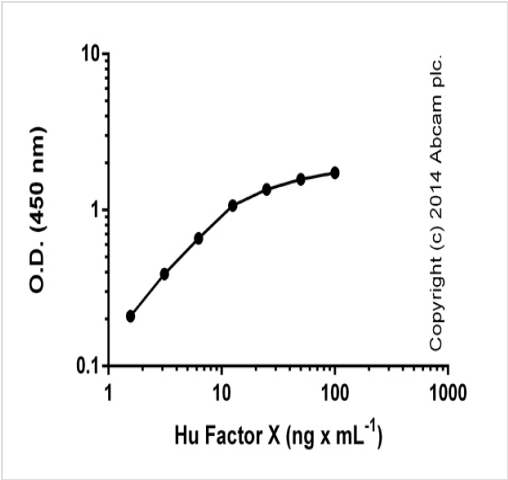
**Tissue specificity** Plasma; synthesized in the liver.

**Involvement in disease** Defects in F10 are the cause of factor X deficiency (FA10D) [MIM:227600]. A hemorrhagic disease with variable presentation. Affected individuals can manifest prolonged nasal and mucosal hemorrhage, menorrhagia, hematuria, and occasionally hemarthrosis. Some patients do not have clinical bleeding diathesis.

**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.

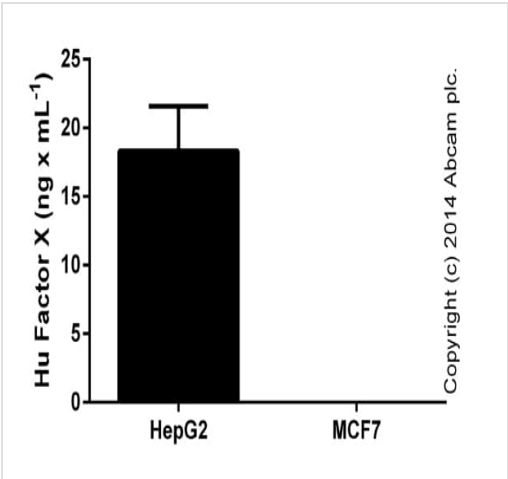
**Post-translational modifications** The vitamin K-dependent, enzymatic carboxylation of some glutamate residues allows the modified protein to bind calcium.  
N- and O-glycosylated.  
The activation peptide is cleaved by factor IXa (in the intrinsic pathway), or by factor VIIa (in the extrinsic pathway).  
The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.

**Cellular localization** Secreted.



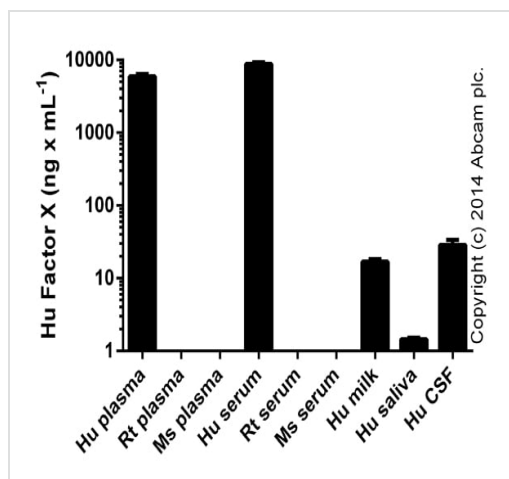
ELISA: Factor X Human ELISA kit (ab108832)

Standard curve with background signal subtracted (duplicates; +/- SD).



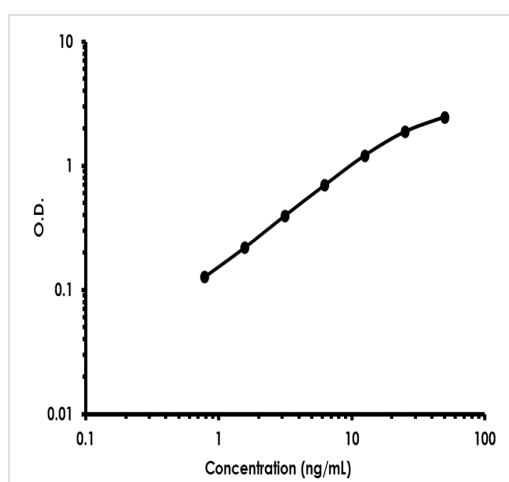
ELISA: Factor X Human ELISA kit (ab108832)

Factor X measured in undiluted culture supernatants with MCF7 levels below level of detection (duplicates +/- SD).



ELISA: Factor X Human ELISA kit (ab108832)

Factor X measured in biological fluids (duplicates +/- SD). Serum and plasma were tested at 1/300-1/3000; other samples at 1/1-1/30.



Typical Standard Curve

Representative Standard Curve using ab108832

**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