

Human Factor X ELISA Kit ab108832

2 References 4 图像

概述

产品名称	人Factor X ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				3.8%
	批次间				
	样品	n	Mean	SD	CV%
	Overall				8.8%
样品类型	Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid				
检测类型	Sandwich (quantitative)				
灵敏度	= 1.5 ng/ml				
范围	0.781 ng/ml - 50 ng/ml				
回收率	97.5 %				
检测时间	4h 00m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's Human Factor <i>Xin vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed				

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.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

功能 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.

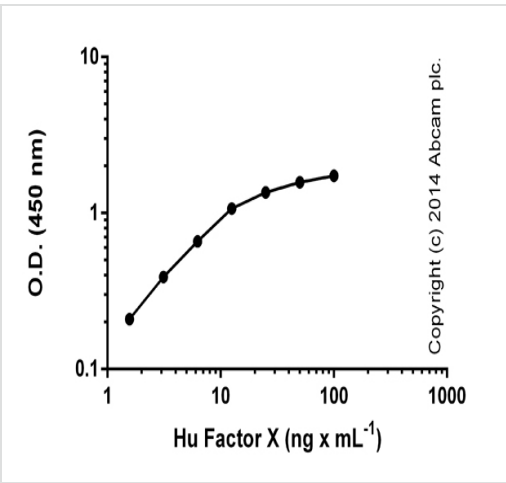
组织特异性 Plasma; synthesized in the liver.

疾病相关 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.

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

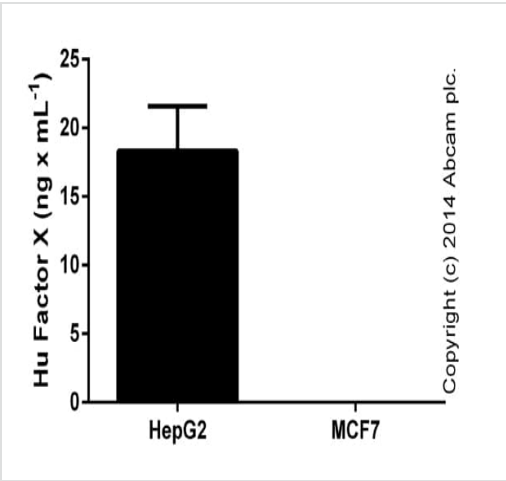
翻译后修饰 The vitamin K-dependent, enzymatic carboxylation of some glutamate residues allows the modified protein to bind calcium.
N- and O-glycosylated.
The activation peptide is cleaved by factor IXa (in the intrinsic pathway), or by factor VIIa (in the extrinsic pathway).
The iron and 2-oxoglutarate dependent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

细胞定位 Secreted.



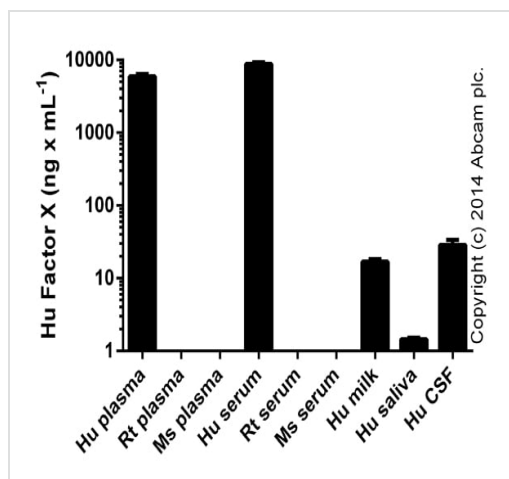
ELISA: Factor X Human ELISA kit (ab108832)

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



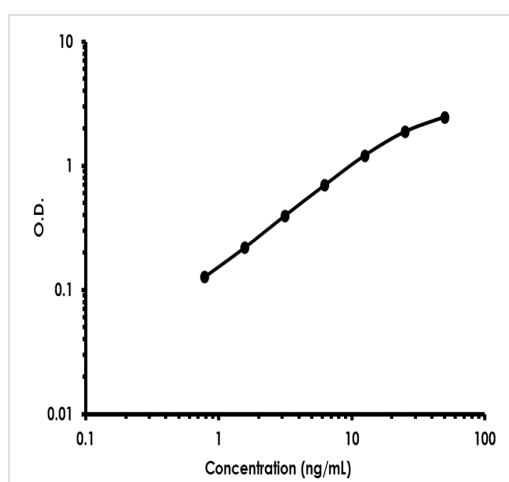
ELISA: Factor X Human ELISA kit (ab108832)

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.



Representative Standard Curve using ab108832

Typical Standard Curve

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors