



Product datasheet

Human Factor XI ELISA Kit ab108834

1 References 1 Image

Overview

Product name	Human Factor XI ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				4.9%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				10%
Sample type	Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	0.69 ng/ml				
Range	1.563 ng/ml - 50 ng/ml				
Recovery	96 %				
Assay time	4h 00m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's Factor XI Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor XI levels in human plasma and serum samples.				

A Factor XI specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor XI 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 XI 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 M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Factor XI Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Factor XI Microplate (12 x 8 well strips)	1 unit
Factor XI Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function

Factor XI triggers the middle phase of the intrinsic pathway of blood coagulation by activating factor IX.

Tissue specificity

Isoform 2 is produced by platelets and megakaryocytes but absent from other blood cells.

Involvement in disease

Defects in F11 are the cause of factor XI deficiency (FA11D) [MIM:612416]; also known as plasma thromboplastin antecedent deficiency or Rosenthal syndrome. It is a hemorrhagic disease characterized by reduced levels and activity of factor XI resulting in moderate bleeding symptoms, usually occurring after trauma or surgery. Patients usually do not present spontaneous bleeding but women can present with menorrhagia. Hemorrhages are usually moderate.

Sequence similarities

Belongs to the peptidase S1 family. Plasma kallikrein subfamily.
Contains 4 apple domains.
Contains 1 peptidase S1 domain.

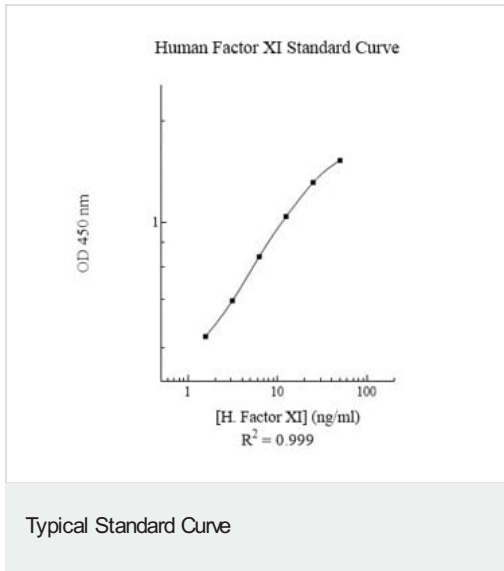
Post-translational modifications

Activated by factor XIIa (or XII), which cleaves each polypeptide after Arg-387 into the light chain, which contains the active site, and the heavy chain, which associates with high molecular weight (HMW) kininogen.

Cellular localization

Secreted.

Images



Representative Standard Curve using ab108834.

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