

Human Factor XIII ELISA Kit ab108836

3 References 画像数 1

製品の概要

製品名	Human Factor XIII ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.4%
サンプルの種類	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 1.5 ng/ml			
検出範囲	5 ng/ml - 80 ng/ml			
添加回収試験	99 %			
全工程の試験時間	4h 0m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Factor XIII Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of factor XIII in plasma, serum, milk, urine, saliva, cell culture, and cell lysate samples.			

A Factor XIII specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor XIII 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 XIII

captured in plate.

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
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 XIII Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Factor XIII Microplate (12 x 8 well strips)	1 unit
Factor XIII Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Factor XIII is activated by thrombin and calcium ion to a transglutaminase that catalyzes the formation of gamma-glutamyl-epsilon-lysine cross-links between fibrin chains, thus stabilizing the fibrin clot. Also cross-link alpha-2-plasmin inhibitor, or fibronectin, to the alpha chains of fibrin.

関連疾患

Defects in F13A1 are the cause of factor XIII subunit A deficiency (FA13AD) [MIM:613225]. FA13AD is an autosomal recessive disorder characterized by a life-long bleeding tendency, impaired wound healing and spontaneous abortion in affected women.

配列類似性

Belongs to the transglutaminase superfamily. Transglutaminase family.

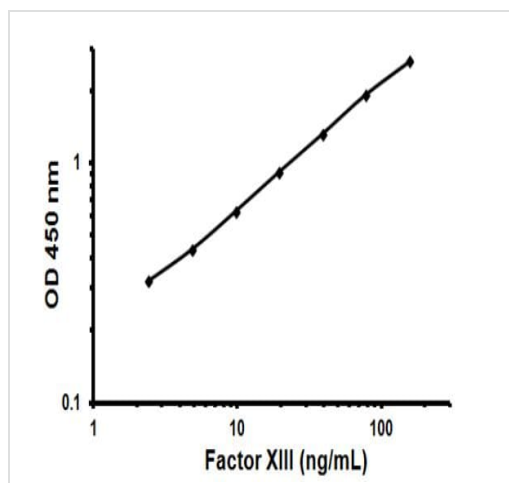
翻訳後修飾

The activation peptide is released by thrombin.

細胞内局在

Cytoplasm. Secreted. Secreted into the blood plasma. Cytoplasmic in most tissues, but also secreted in the blood plasma.

画像



Representative Standard Curve using ab108836

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.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