

Human Factor V ELISA Kit ab137976

2 References 画像数 1

製品の概要

製品名	Human Factor V 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	CV%
	Overall				4.8%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Overall				10.9%
サンプルの種類	Cell culture supernatant, Milk, Urine, Plasma, Cerebral Spinal Fluid				
アッセイタイプ	Sandwich (quantitative)				
検出感度	> 49 pg/ml				
検出範囲	2.5 ng/ml - 10 ng/ml				
添加回収試験	102 %				
	特定サンプルでの回収試験				
	サンプルの種類	平均 %		測定範囲	
	Milk			91% - 103%	
	Urine			94% - 102%	
	Plasma			94% - 104%	
全工程の試験時間	4h 00m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's Factor V Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor V in plasma, urine, milk, and cell culture supernatants.				

A Factor V specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor V specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate 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 V captured in plate.

Factor V Standard is a native protein.

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

機能

Central regulator of hemostasis. It serves as a critical cofactor for the prothrombinase activity of factor Xa that results in the activation of prothrombin to thrombin.

組織特異性

Plasma.

関連疾患

Factor V deficiency
Thrombophilia due to activated protein C resistance
Budd-Chiari syndrome
Ischemic stroke
Pregnancy loss, recurrent, 1

配列類似性

Belongs to the multicopper oxidase family.

Contains 3 F5/8 type A domains.
Contains 2 F5/8 type C domains.
Contains 6 plastocyanin-like domains.

ドメイン

Domain B contains 35 x 9 AA tandem repeats, and 2 x 17 AA repeats.

翻訳後修飾

Thrombin activates factor V proteolytically to the active cofactor, factor Va (formation of a heavy chain at the N-terminus and a light chain at the C-terminus).

Sulfation is required for efficient thrombin cleavage and activation and for full procoagulant activity.

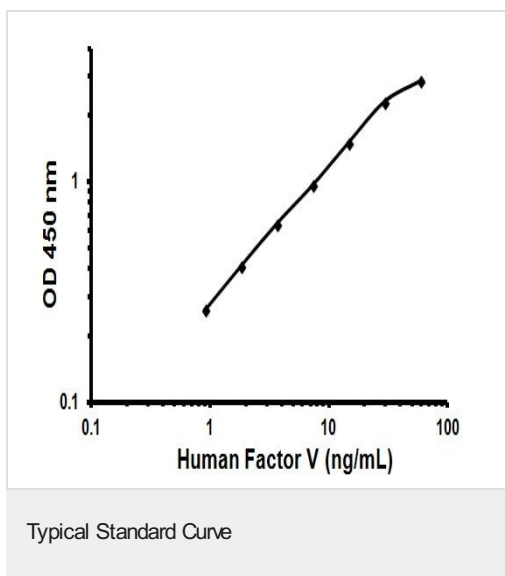
Activated protein C inactivates factor V and factor Va by proteolytic degradation.

Phosphorylation sites are present in the extracellular medium.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab137976

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