

### Human Factor VII ELISA Kit ab108829

5 References [画像数 1](#)

#### 製品の概要

<b>製品名</b>	Human Factor VII ELISA Kit										
<b>検出方法</b>	Colorimetric										
<b>再現性</b>	Intra-Assay (同時再現性)										
	<table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="width: 25%;">サンプル</th> <th style="width: 10%;">N</th> <th style="width: 25%;">平均値</th> <th style="width: 15%;">SD</th> <th style="width: 25%;">CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>4.9%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				4.9%
サンプル	N	平均値	SD	CV%							
Overall				4.9%							
	Inter-Assay (日差再現性)										
	<table border="1" style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="width: 25%;">サンプル</th> <th style="width: 10%;">N</th> <th style="width: 25%;">平均値</th> <th style="width: 15%;">SD</th> <th style="width: 25%;">CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>8.6%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				8.6%
サンプル	N	平均値	SD	CV%							
Overall				8.6%							
<b>サンプルの種類</b>	Cell culture supernatant, Saliva, Serum, Plasma										
<b>アッセイタイプ</b>	Sandwich (quantitative)										
<b>検出感度</b>	= 1.8 ng/ml										
<b>検出範囲</b>	10 ng/ml - 40 ng/ml										
<b>添加回収試験</b>	96 %										
<b>全工程の試験時間</b>	4h 00m										
<b>ステップ</b>	Multiple steps standard assay										
<b>種交差性</b>	交差種: Human										
<b>製品の概要</b>	<p>Abcam's Factor VII Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor VII concentrations in Human cell culture supernatants, serum and plasma.</p> <p>A Factor VII specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor VII 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 VII captured in plate. This assay recognizes FVII and FVIIa.</p>										

Get higher sensitivity in only 90 minutes with Human Factor VII ELISA Kit (**ab190810**) from our SimpleStep ELISA® range.

**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 N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Factor VII Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Factor VII Microplate (12 x 8 well strips)	1 unit
Factor VII Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Initiates the extrinsic pathway of blood coagulation. Serine protease that circulates in the blood in a zymogen form. Factor VII is converted to factor VIIa by factor Xa, factor XIIa, factor IXa, or thrombin by minor proteolysis. In the presence of tissue factor and calcium ions, factor VIIa then converts factor X to factor Xa by limited proteolysis. Factor VIIa will also convert factor IX to factor IXa in the presence of tissue factor and calcium.

組織特異性

Plasma.

関連疾患

Defects in F7 are the cause of factor VII deficiency (FA7D) [MIM:227500]. FA7D is a rare hereditary hemorrhagic disease. The clinical picture can be very severe, with the early occurrence of intracerebral hemorrhages or hemarthroses, or, in contrast, moderate with cutaneous-mucosal hemorrhages (epistaxis, menorrhagia) or hemorrhages provoked by a surgical intervention. Numerous subjects are completely asymptomatic despite a very low F7 level.

配列類似性

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

翻訳後修飾

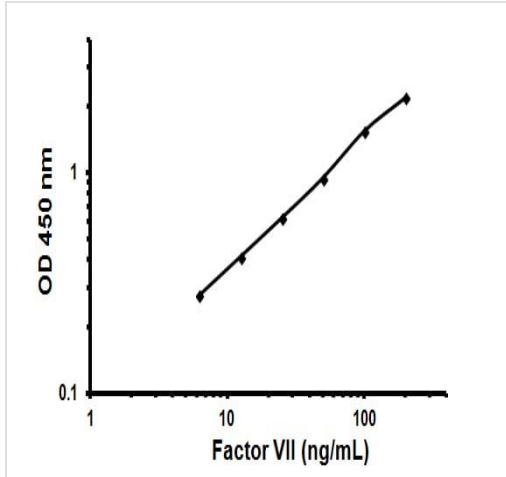
The vitamin K-dependent, enzymatic carboxylation of some glutamate residues allows the modified protein to bind calcium.

The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.

細胞内局在

Secreted.

## 画像



Representative Standard Curve using ab108829.

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