

Human Factor VII ELISA Kit ab108829

5 References 1 图像

概述

产品名称	人Factor VII ELISA试剂盒																				
检测方法	Colorimetric																				
精确度	<div>批次内</div> <table><tr><td>样品</td><td>n</td><td>Mean</td><td>SD</td><td>CV%</td></tr><tr><td>Overall</td><td></td><td></td><td></td><td>4.9%</td></tr></table> <div>批次间</div> <table><tr><td>样品</td><td>n</td><td>Mean</td><td>SD</td><td>CV%</td></tr><tr><td>Overall</td><td></td><td></td><td></td><td>8.6%</td></tr></table>	样品	n	Mean	SD	CV%	Overall				4.9%	样品	n	Mean	SD	CV%	Overall				8.6%
样品	n	Mean	SD	CV%																	
Overall				4.9%																	
样品	n	Mean	SD	CV%																	
Overall				8.6%																	
样品类型	Cell culture supernatant, Saliva, Serum, Plasma																				
检测类型	Sandwich (quantitative)																				
灵敏度	= 1.8 ng/ml																				
范围	10 ng/ml - 40 ng/ml																				
回收率	96 %																				
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 Factor VII Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed																				

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.

平台 Microplate

性能

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

组件	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

功能 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.

组织特异性 Plasma.

疾病相关 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.

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.

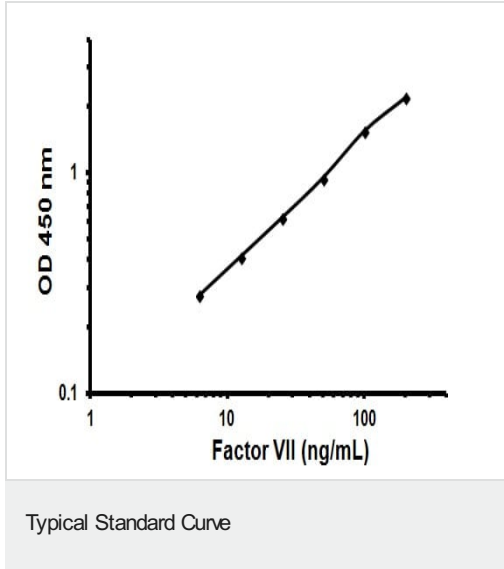
翻译后修饰 The vitamin K-dependent, enzymatic carboxylation of some glutamate residues allows the modified protein to bind calcium.

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

细胞定位

Secreted.

图片



Representative Standard Curve using ab108829.

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