

# Factor VII Human Chromogenic Activity Assay Kit ab108830

[1 References](#) [2 Images](#)

Overview

Product name	Factor VII Human Chromogenic Activity Assay Kit																							
Detection method	Colorimetric																							
Precision	<div>Intra-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td>60</td><td></td><td></td><td>5.2%</td></tr></table> <div>Inter-assay</div> <table><tr><th>Sample</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td>60</td><td></td><td></td><td>10.5%</td></tr></table>				Sample	n	Mean	SD	CV%	Overall	60			5.2%	Sample	n	Mean	SD	CV%	Overall	60			10.5%
Sample	n	Mean	SD	CV%																				
Overall	60			5.2%																				
Sample	n	Mean	SD	CV%																				
Overall	60			10.5%																				
Sample type	Cell culture supernatant, Serum, Plasma																							
Assay type	Enzyme activity (quantitative)																							
Sensitivity	5 mIU/ml																							
Assay time	3h 00m																							
Species reactivity	<b>Reacts with:</b> Human																							
Product overview	<p>Factor VII Human Chromogenic Activity Assay Kit (ab108830) has been developed to determine human FVII activity in plasma, serum and cell culture supernatants. The assay couples immunofunctional and indirect amidolytic assay. A polyclonal antibody specific for Human FVII has been pre-coated onto a microplate and active FVII is bound to the immobilized antibody. The assay measures the ability of lipoprotein TF/FVIIa to activate factor X (FX) to factor Xa. The amidolytic activity of the TF/FVIIa complex is quantitated by the amount of FXa produced using a highly specific FXa substrate releasing a yellow para-nitroaniline (pNA) chromophore. The change in absorbance of the pNA at 405 nm is directly proportional to the FVII enzymatic activity.</p> <p><b>The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.</b></p>																							

Notes	Factor VII (FVII) is a vitamin K-dependent plasma glycoprotein that is synthesized in the liver and circulates in blood as a single-chain inactive zymogen with a molecular mass of 50 kDa. Upon tissue damage and vascular injury, the cell surface receptor and cofactor tissue factor (TF) binds
-------	---

and allosterically activates FVII to its active form, FVIIa. The TF/FVIIa complex catalyzes the conversion of both factor IX to factor IXa and factor X to factor Xa to initiate coagulation via the extrinsic pathway. Very low levels of FVII are associated with severe coagulation disorders. Elevated plasma levels of FVII coagulant activity constitute an independent risk factor for fatal outcomes of coronary heart disease in middle-aged men.

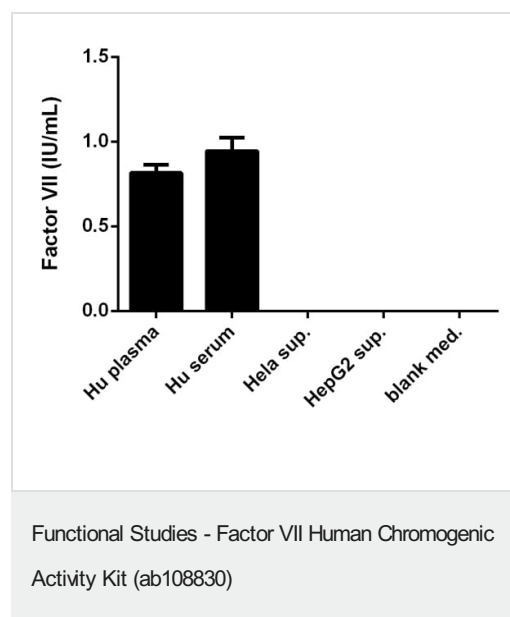
**Platform** Microplate reader

## Properties

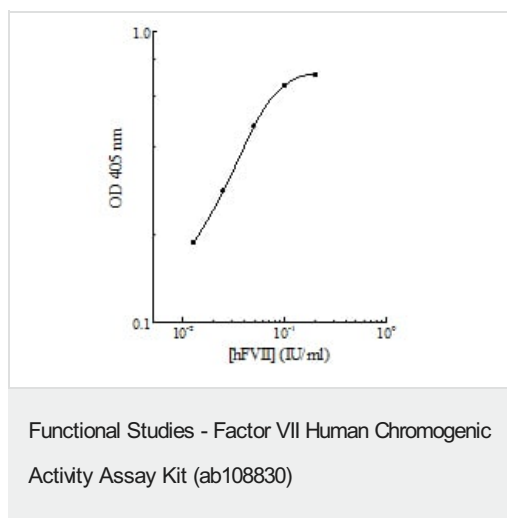
**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
10X Diluent M Concentrate	1 x 20ml
20X Wash Buffer Concentrate	1 x 30ml
Assay Diluent	1 x 20ml
FXa Substrate	2 units
Human FVII Microplate	1 x 96 units
Human FVII Standard	1 unit
Human FX	1 unit
rhTF (lipoprotein)	1 unit
Sealing Tapes	3 units

## Images



Factor VII measured in biological fluids showing quantity (IU) per mL of tested sample



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.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