

Product datasheet

Factor VIIIa Activity Assay Kit (Fluorometric) ab204696

[2 Images](#)

Overview

Product name	Factor VIIIa Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Serum, Plasma, Purified protein
Assay type	Enzyme activity
Sensitivity	1 ng/well
Product overview	<p>Factor VIIIa Activity Assay Kit (Fluorometric) (ab204696) is based on the ability of FVIIIa to generate FXa. The generated FXa proteolytically cleaves a synthetic substrate and releases a fluorophore, AMC, which can be easily quantified by fluorescence microplate reader.</p> <p>The assay is simple, rapid and can detect activity as low as 1 ng of FVIIIa in a variety of samples.</p>
Notes	<p>The coagulation Factor VIII (anti-hemophilic factor, AHF) is a vital blood-clotting protein. Factor VIII circulates in the bloodstream as an inactive protein, bound to a large multimeric glycoprotein called von Willebrand factor (VWF). Upon an injury to the blood vessel, VWF dissociates from FVIII and releases the active form of FVIII (FVIIIa). In the presence of calcium ions and negatively charged membrane phospholipids, activated factor VIII (FVIIIa) then binds to the activated Factor IX (FIXa) and proteolytically activates factor X (FX) to factor Xa (FXa).</p>
Platform	Microplate reader

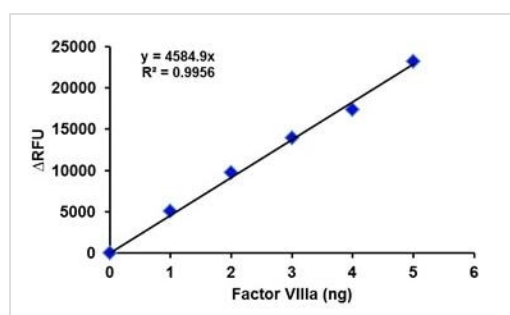
Properties

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

Components	100 tests
Enzyme Mix I	1 vial
Enzyme Mix II	1 vial
FVIIIa Assay Buffer	1 x 15ml
FVIIIa Enzyme Standard	1 x 2.6µg
FXa Substrate-AMC	1 x 200µl

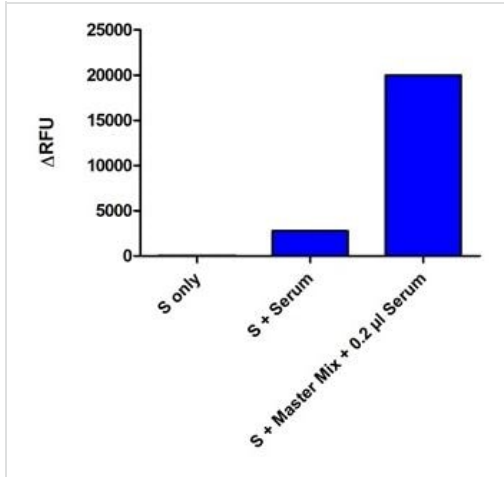
Components	100 tests
Phospholipids	1 x 600µl
Function	Factor VIII, along with calcium and phospholipid, acts as a cofactor for factor IXa when it converts factor X to the activated form, factor Xa.
Involvement in disease	Defects in F8 are the cause of hemophilia A (HEMA) [MIM:306700]. A disorder of blood coagulation characterized by a permanent tendency to hemorrhage. About 50% of patients have severe hemophilia resulting in frequent spontaneous bleeding into joints, muscles and internal organs. Less severe forms are characterized by bleeding after trauma or surgery. Note=Of particular interest for the understanding of the function of F8 is the category of CRM (cross-reacting material) positive patients (approximately 5%) that have considerable amount of F8 in their plasma (at least 30% of normal), but the protein is non-functional; i.e., the F8 activity is much less than the plasma protein level. CRM-reduced is another category of patients in which the F8C antigen and activity are reduced to approximately the same level. Most mutations are CRM negative, and probably affect the folding and stability of the protein.
Sequence similarities	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.
Domain	Domain F5/8 type C 2 is responsible for phospholipid-binding and essential for factor VIII activity.
Post-translational modifications	Sulfation on Tyr-1699 is essential for binding vWF.
Cellular localization	Secreted > extracellular space.

Images



Typical Factor VIIIa Standard Curve using Factor VIIIa Activity Assay Kit (Fluorometric) (ab204696).

Factor VIIIa Activity Assay Kit (Fluorometric)
(ab204696)



Factor VIIIa activity was measured in serum samples in the presence and absence of the master mix. S: Substrate.

Factor VIIIa Activity Assay Kit (Fluorometric)
(ab204696)

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