

Factor IXa Activity Assay Kit (Fluorometric) ab204727

[1 References](#) [2 图像](#)

概述

产品名称	Factor IXa Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Serum, Plasma, Purified protein
检测类型	Enzyme activity
产品概述	Factor IXa Activity Assay Kit (Fluorometric) (ab204727) is based on the ability of FIXa 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. The assay is simple, rapid and can detect activity as low as 10 pg of FIXa in a variety of samples.

说明

This product is manufactured by BioVision, an Abcam company and was previously called K364 Factor IXa Activity Assay Kit (Fluorometric). K364-100 is the same size as the 100 test size of ab204727.

The coagulation Factor IX (or Christmas factor, EC 3.4.21.22) is a vitamin K-dependent serine protease. Factor IX is produced as an inactive precursor and is activated via cleavage by either factor XIa (contact pathway) or factor VIIa (tissue factor pathway). In the presence of calcium ions and negatively charged membrane phospholipids, activated factor IX (FIXa) then binds to the activated Factor VIII (FVIIIa) and proteolytically activates factor X (FX) to factor Xa (FXa).

平台 Microplate reader

性能

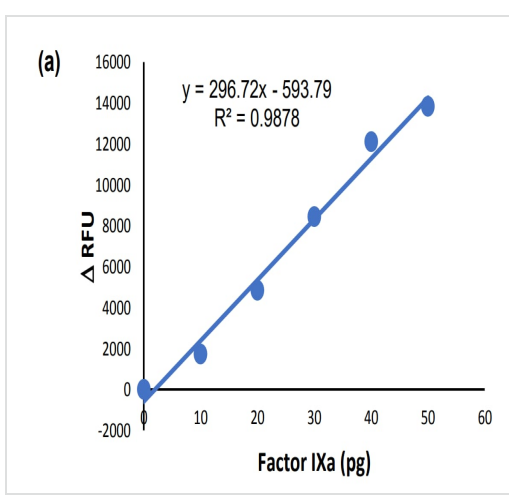
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
Active Factor VIIIa	1 vial
Assay Buffer XLIV	1 x 15ml
Enzyme Mix XXII	1 vial
FIXa Enzyme Standard	1 x 10ng
FXa Substrate	1 x 200µl

组件	100 tests
Phospholipid Mixture	1 x 600µl

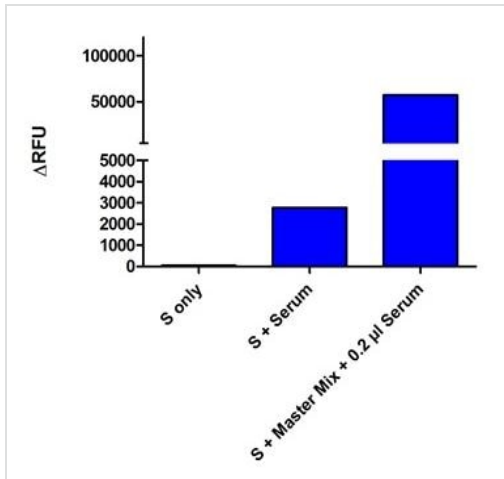
功能	Factor IX is a vitamin K-dependent plasma protein that participates in the intrinsic pathway of blood coagulation by converting factor X to its active form in the presence of Ca(2+) ions, phospholipids, and factor VIIIa.
组织特异性	Synthesized primarily in the liver and secreted in plasma.
疾病相关	Defects in F9 are the cause of recessive X-linked hemophilia B (HEMB) [MIM:306900]; also known as Christmas disease. Note=Mutations in position 43 (Oxford-3, San Dimas) and 46 (Cambridge) prevents cleavage of the propeptide, mutation in position 93 (Alabama) probably fails to bind to cell membranes, mutation in position 191 (Chapel-Hill) or in position 226 (Nagoya OR Hilo) prevent cleavage of the activation peptide. Defects in F9 are the cause of thrombophilia due to factor IX defect (THR-FIX) [MIM:300807]. A hemostatic disorder characterized by a tendency to thrombosis.
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.
结构域	Calcium binds to the gamma-carboxyglutamic acid (Gla) residues and, with stronger affinity, to another site, beyond the Gla domain.
翻译后修饰	Activated by factor XIa, which excises the activation peptide. 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



Typical Factor IXa Standard Curve.

Factor IX Activity Assay Kit (Fluorometric)
(ab204727)



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

Factor IX Activity Assay Kit (Fluorometric)
(ab204727)

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