

## 1 References 画像数 2

**医藥用外劇物**

製品名	Factor IXa Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Serum, Plasma, Purified protein
アッセイタイプ	Enzyme activity
製品の概要	<p>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.</p>
特記事項	<p>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.</p> <p>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).</p>
試験プラットフォーム	Microplate reader

## 製品の特長

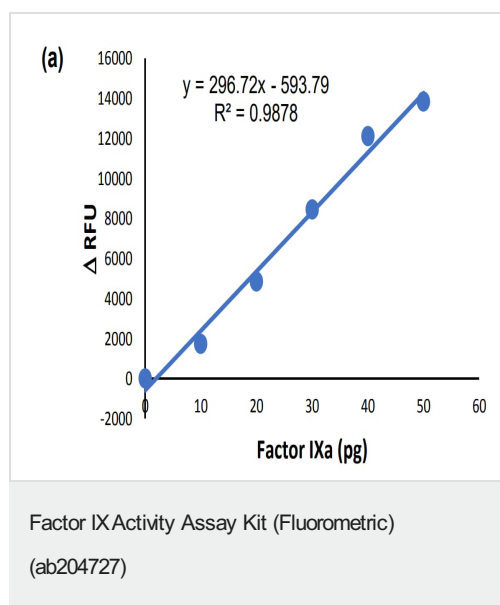
**保存方法** Store at -20°C. Please refer to protocols.

内容	100 tests
Active Factor VIIIa	1 vial
Assay Buffer XLIV	1 x 15ml
Enzyme Mix XXII	1 vial
FIXa Enzyme Standard	1 x 10ng

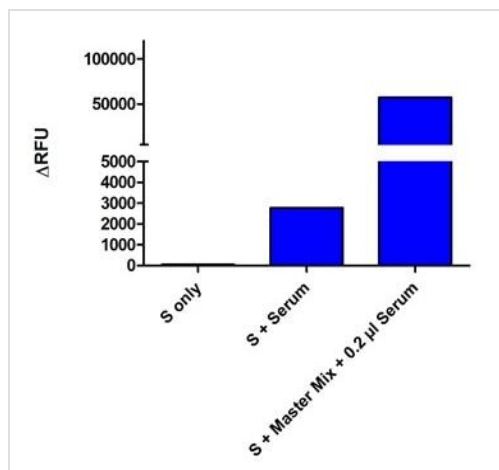
内容	100 tests
FXa Substrate	1 x 200μl
Phospholipid Mixture	1 x 600μl

機能	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.
組織特異性	Synthesized primarily in the liver and secreted in plasma.
関連疾患	<p>Defects in F9 are the cause of recessive X-linked hemophilia B (HEMB) [MIM:306900]; also known as Christmas disease.</p> <p>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.</p> <p>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.</p>
配列類似性	<p>Belongs to the peptidase S1 family.</p> <p>Contains 2 EGF-like domains.</p> <p>Contains 1 Gla (gamma-carboxy-glutamate) domain.</p> <p>Contains 1 peptidase S1 domain.</p>
ドメイン	Calcium binds to the gamma-carboxyglutamic acid (Gla) residues and, with stronger affinity, to another site, beyond the Gla domain.
翻訳後修飾	<p>Activated by factor XIa, which excises the activation peptide.</p> <p>The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.</p>
細胞内局在	Secreted.

## 画像



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

**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