

Factor Xla Assay Kit ab241037

4 图像

概述

产品名称	Factor Xla Assay试剂盒
检测方法	Colorimetric
产品概述	<p>The Factor Xla Assay Kit (ab241037) utilizes the ability of factor Xla to cleave a synthetic substrate to release <i>p</i>-Nitroaniline (<i>p</i>NA) which can be quantitatively measured by a colorimetric assay (OD 405 nm).</p> <p>The kit is easy-to-use and can detect Factor Xla (as low as 1 mPEU) from plasma and purified protein samples.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K973 Factor Xla Activity Assay Kit (Colorimetric). K973-100 is the same size as the 100 test size of ab241037.</p>

性能

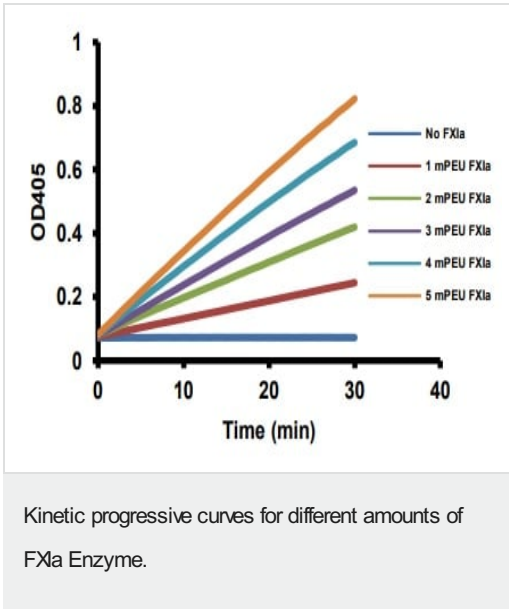
存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	100 tests
Assay Buffer XXXV	1 x 25ml
FXla Activator	1 x 1ml
FXla Substrate	1 x 1ml
FXIIa Inhibitor	1 x 0.1ml
Human Factor Xla	1 vial
PK Inhibitor	1 x 0.1ml
pNA Standard I	1 x 20µl

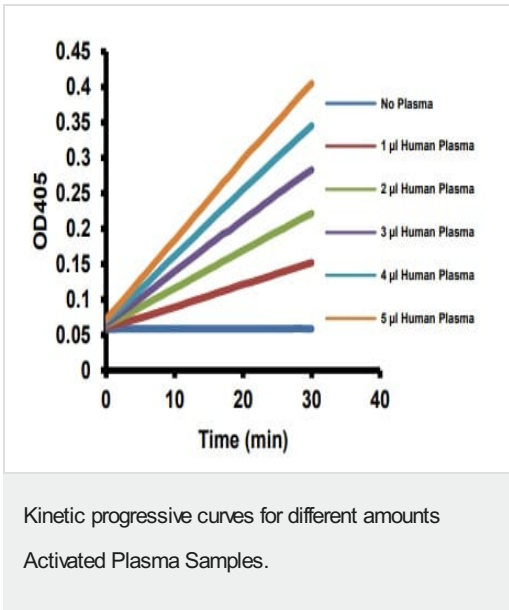
功能	Factor XI triggers the middle phase of the intrinsic pathway of blood coagulation by activating factor IX.
----	--

组织特异性	Isoform 2 is produced by platelets and megakaryocytes but absent from other blood cells.
疾病相关	Defects in F11 are the cause of factor XI deficiency (FA11D) [MIM:612416]; also known as plasma thromboplastin antecedent deficiency or Rosenthal syndrome. It is a hemorrhagic disease characterized by reduced levels and activity of factor XI resulting in moderate bleeding symptoms, usually occurring after trauma or surgery. Patients usually do not present spontaneous bleeding but women can present with menorrhagia. Hemorrhages are usually moderate.
序列相似性	Belongs to the peptidase S1 family. Plasma kallikrein subfamily. Contains 4 apple domains. Contains 1 peptidase S1 domain.
翻译后修饰	Activated by factor XIIa (or XII), which cleaves each polypeptide after Arg-387 into the light chain, which contains the active site, and the heavy chain, which associates with high molecular weight (HMW) kininogen.
细胞定位	Secreted.

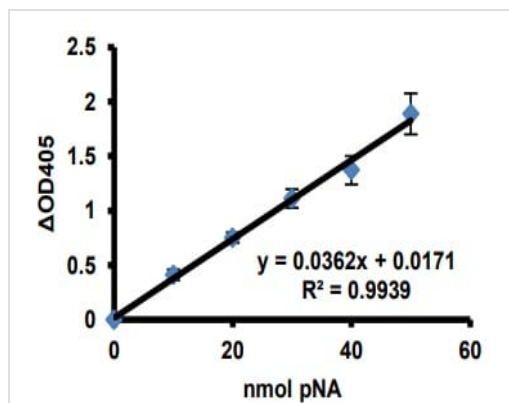
图片



Kinetic progressive curves for different amounts of FXIa Enzyme.

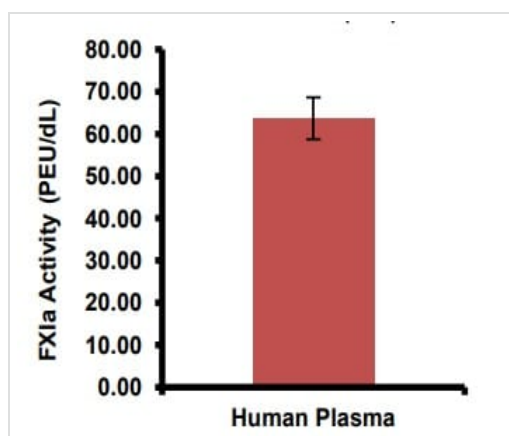


Kinetic progressive curves for different amounts Activated Plasma Samples.



Standard curve for pNA.

Standard curve for pNA.



Standard curve for pNA (n = 3) was used to estimate FXIa activity in plasma (n = 5). Assays were performed according to the kit protocol.

FXIa activity in plasma.

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