

Factor Xa Inhibitor Screening Assay Kit (Fluorometric)  
ab204712

2 Images

Overview

Product name	Factor Xa Inhibitor Screening Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Inhibitor compounds
Assay type	Enzyme activity
Product overview	Factor Xa Inhibitor Screening Assay Kit (Fluorometric) (ab204712) is based on the ability of Factor Xa to cleave a synthetic substrate thereby releasing a fluorophore, AMC, which can be quantified by a fluorescence reader at Ex/Em = 350/450 nm. In the presence of an inhibitor, the extent of the cleavage reaction is reduced or completely abolished. The loss in the fluorescence intensity can be correlated to the amount of inhibitor present in the assay solution. The Factor Xa Inhibitor Screening Kit is a simple, straightforward, high-throughput assay to screen Factor Xa inhibitors.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K362 Factor Xa Inhibitor Screening Kit (Fluorometric). K362-100 is the same size as the 100 test size of ab204712.</p> <p>Factor Xa (FXa) is the activated form of the coagulation factor X (Stuart-Power factor, thrombokinase, prothrombinase, thromboplastin, E.C.3.4.21.6). Factor X, a serine endopeptidase, plays an important role at several stages of the coagulation pathway. It acts by converting prothrombin into active thrombin by complexing with activated co-factor V in the prothrombinase complex. Unfractionated heparin and various low molecular weight heparins bind to plasma cofactor Antithrombin to inactivate several coagulation factors including factor Xa.</p>
Platform	Microplate reader

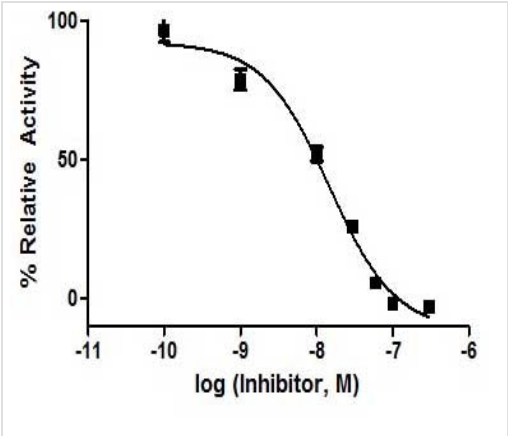
Properties

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

Components	100 tests
FXa Assay Buffer	1 x 15ml
FXa Dilution Buffer	1 x 1ml

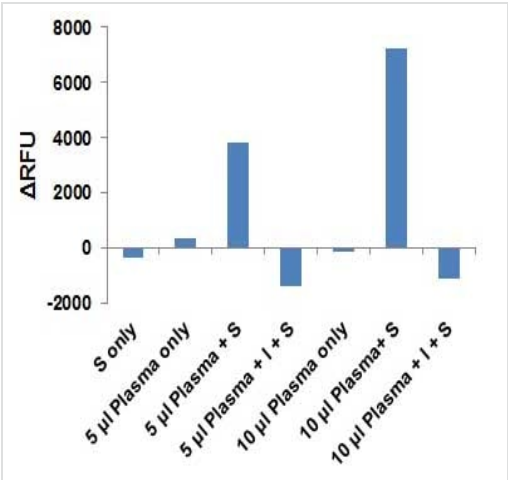
Components	100 tests
FXa Enzyme	1 x 5µl
FXa Inhibitor	1 x 10µl
FXa Substrate	1 x 200µl

Images



Inhibition of Factor Xa activity by FXa Inhibitor (GGACK Dihydrochloride).

Factor Xa Inhibitor Screening Assay Kit  
(Fluorometric) (ab204712)



Factor Xa activity was measured in plasma samples in the presence and absence of FXa Inhibitor (GGACK Dihydrochloride).

S = Substrate, I = Inhibitor.

Factor Xa Inhibitor Screening Assay Kit  
(Fluorometric) (ab204712)

**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