

Thrombin Activity Assay Kit (Fluorometric) ab197006

5 References [画像数 2](#)

製品の概要

製品名	Thrombin Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Plasma, Purified protein
アッセイタイプ	Quantitative
検出感度	1 ng
種交差性	交差種: Mammals, Other species
製品の概要	Thrombin Activity Assay Kit (Fluorometric) (ab197006) is simple, rapid and can detect Thrombin activity as low as 1 ng in samples.

In the thrombin activity assay protocol, a synthetic AMC-based peptide substrate is proteolytically cleaved by thrombin to release a fluorophore, AMC, which can be easily quantified by fluorescence reader at Ex/Em = 350/450 nm.

Thrombin activity assay protocol summary:
 - add standards and samples to wells
 - add reaction mix
 - analyze with a microplate reader for 30-60 min

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K373 Thrombin Activity Fluorometric Assay Kit. K373-100 is the same size as the 100 test size of ab197006.</p> <p>Thrombin enzyme (Factor IIa, EC 3.4.21.5), a serine protease, is an important clotting factor in the coagulation cascade that involves the conversion of soluble fibrinogen to insoluble active fibrin strands.</p>
------	--

試験プラットフォーム	Microplate reader
------------	-------------------

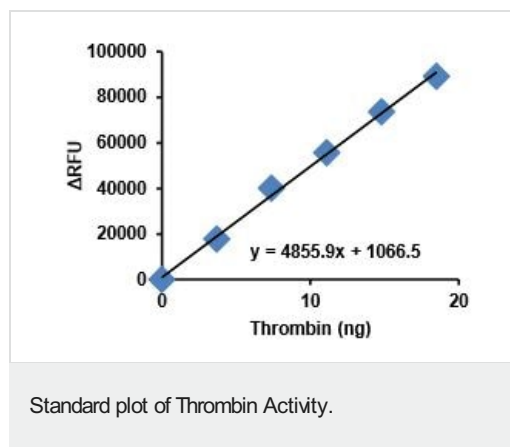
製品の特性

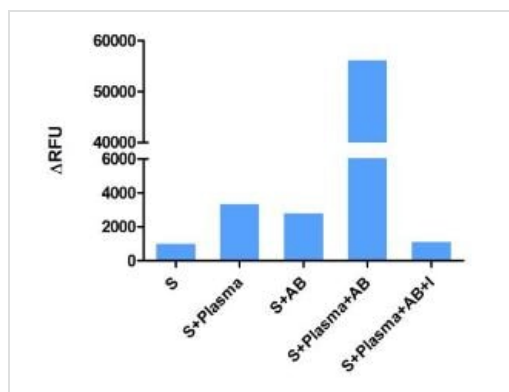
保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	100 tests
Thrombin Assay Buffer	1 x 15ml
Thrombin Dilution Buffer	1 x 1ml
Thrombin Standard	1 x 5μl
Thrombin Substrate	1 x 0.5ml

機能	Thrombin, which cleaves bonds after Arg and Lys, converts fibrinogen to fibrin and activates factors V, VII, VIII, XIII, and, in complex with thrombomodulin, protein C. Functions in blood homeostasis, inflammation and wound healing.
組織特異性	Expressed by the liver and secreted in plasma.
関連疾患	Factor II deficiency Ischemic stroke Thrombophilia due to thrombin defect Pregnancy loss, recurrent, 2
配列類似性	Belongs to the peptidase S1 family. Contains 1 Gla (gamma-carboxy-glutamate) domain. Contains 2 kringle domains. Contains 1 peptidase S1 domain.
翻訳後修飾	The gamma-carboxyglutamyl residues, which bind calcium ions, result from the carboxylation of glutamyl residues by a microsomal enzyme, the vitamin K-dependent carboxylase. The modified residues are necessary for the calcium-dependent interaction with a negatively charged phospholipid surface, which is essential for the conversion of prothrombin to thrombin. N-glycosylated. N-glycan heterogeneity at Asn-121: Hex3HexNAc3 (minor), Hex4HexNAc3 (minor) and Hex5HexNAc4 (major). At Asn-143: Hex4HexNAc3 (minor) and Hex5HexNAc4 (major).
細胞内局在	Secreted, extracellular space.

画像





Thrombin Activity in Plasma.

Thrombin activity was measured in plasma samples in the presence and absence of a Thrombin inhibitor, PPACK Dihydrochloride. S = Substrate, I = Inhibitor, AB = Activation Buffer containing Factor Xa. Assays were performed following the kit protocol.

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