

Product datasheet

Plasmin Activity Assay Kit (Fluorometric) ab204728

★★★★★ [1 Abreviews](#) [2 References](#) [2 Images](#)

Overview

Product name	Plasmin Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Plasma, Purified protein
Assay type	Enzyme activity
Sensitivity	10 ng
Assay time	0h 20m
Species reactivity	Reacts with: Mammals, Other species
Product overview	Plasmin Activity Assay Kit (Fluorometric) (ab204728) is based on the ability of Plasmin to proteolytically cleave a synthetic plasmin substrate and release a fluorophore, AMC, which can be easily quantified by fluorescence microplate readers at Ex/Em = 360/450 nm. This assay kit is simple, rapid and can detect plasmin activity as low as 10 ng in a variety of samples.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K381 Plasmin Activity Assay Kit (Fluorometric). K381-100 is the same size as the 100 test size of ab204728.</p> <p>Plasmin (EC 3.4.21.7) is a serine protease occurring in plasma as plasminogen. Upon activation via cleavage by plasminogen activators; plasmin solubilizes fibrin clots and activates and/or degrades compounds of the coagulation and complement systems. Plasminogen-Plasmin system has also been implicated in a wide variety of physiologic and pathologic processes, including tumor growth, invasion and metastasis.</p>
Platform	Microplate reader

Properties

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

Components	100 tests
Plasmin Assay Buffer	1 x 15ml
Plasmin Dilution Buffer	1 x 1.5ml
Plasmin Enzyme	1 x 5µl

Components	100 tests
Plasmin Substrate I	1 x 200µl

Function

Plasmin dissolves the fibrin of blood clots and acts as a proteolytic factor in a variety of other processes including embryonic development, tissue remodeling, tumor invasion, and inflammation. In ovulation, weakens the walls of the Graafian follicle. It activates the urokinase-type plasminogen activator, collagenases and several complement zymogens, such as C1 and C5. Cleavage of fibronectin and laminin leads to cell detachment and apoptosis. Also cleaves fibrin, thrombospondin and von Willebrand factor. Its role in tissue remodeling and tumor invasion may be modulated by CSPG4. Binds to cells.

Angiostatin is an angiogenesis inhibitor that blocks neovascularization and growth of experimental primary and metastatic tumors in vivo.

Tissue specificity

Present in plasma and many other extracellular fluids. It is synthesized in the liver.

Involvement in disease

Plasminogen deficiency

Sequence similarities

Belongs to the peptidase S1 family. Plasminogen subfamily.

Contains 5 kringle domains.

Contains 1 PAN domain.

Contains 1 peptidase S1 domain.

Domain

Kringle domains mediate interaction with CSPG4.

Post-translational modifications

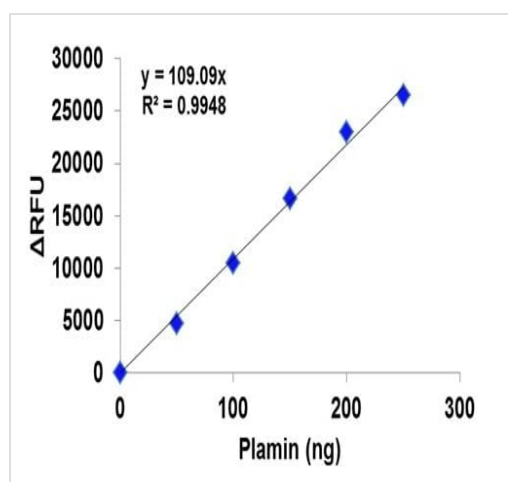
N-linked glycan contains N-acetylglucosamine and sialic acid. O-linked glycans consist of Gal-GalNAc disaccharide modified with up to 2 sialic acid residues (microheterogeneity).

In the presence of the inhibitor, the activation involves only cleavage after Arg-580, yielding two chains held together by two disulfide bonds. In the absence of the inhibitor, the activation involves additionally the removal of the activation peptide.

Cellular localization

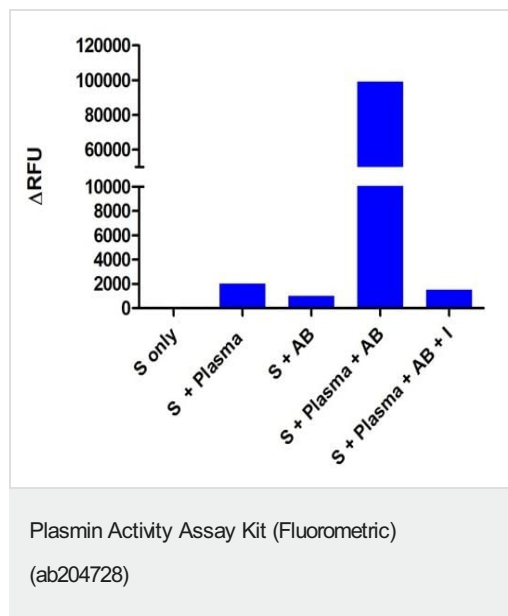
Secreted. Locates to the cell surface where it is proteolytically cleaved to produce the active plasmin. Interaction with HRG tethers it to the cell surface.

Images



Typical Plasmin standard curve.

Plasmin Activity Assay Kit (Fluorometric)
(ab204728)



Plasmin activity was measured in plasma samples in the presence and absence of a Plasmin inhibitor, Aprotinin. S = Substrate, I = Inhibitor, AB = Activation Buffer containing Urokinase

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