

Product datasheet

Plasmin Activity Assay Kit (Colorimetric) ab273301

[3 Images](#)

Overview

Product name	Plasmin Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Plasma
Assay type	Quantitative
Assay duration	Multiple steps standard assay
Product overview	Plasmin Activity Assay Kit (Colorimetric) (ab273301) utilizes the ability of active plasmin to hydrolyze the synthetic substrate thereby releasing pNA (chromophore), which can be easily quantified at OD 405 nm. The stable colorimetric signal is directly proportional to the plasmin activity in samples. The kit includes a specific inhibitor that can be used to compensate for the potential non-specific background in unknown samples. Our assay kit is simple, specific and can detect as low as 40 μ U of plasmin activity in samples.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called K945 Plasmin Activity Assay Kit (Colorimetric). K945-100 is the same size as the 100 test size of ab273301.
Platform	Microplate

Properties

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

Components	100 tests
pNA Standard I	1 x 20 μ l
Assay Buffer A1	1 x 50ml
Plasmin Inhibitor Mix	1 vial
Plasmin Positive Control	1 x 15 μ l
Plasmin Substrate II	1 x 400 μ l

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

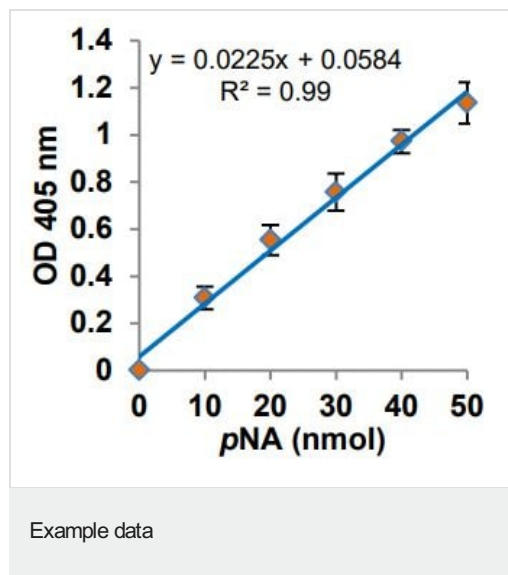
N-linked glycan contains N-acetylglucosamine and sialic acid. O-linked glycans consist of Gal-GalNAc disaccharide modified with up to 2 sialic acid residues (microheterogeneity).

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.

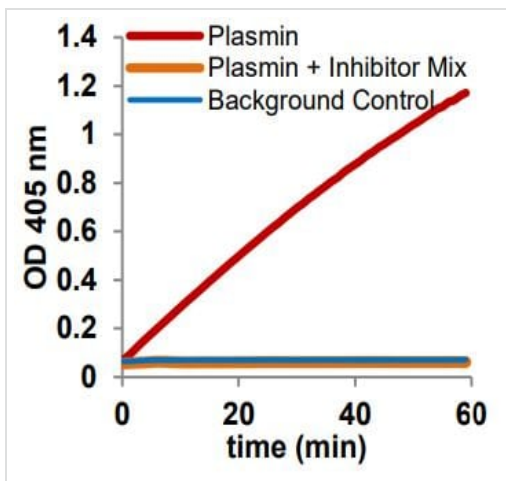
Cellular localization

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

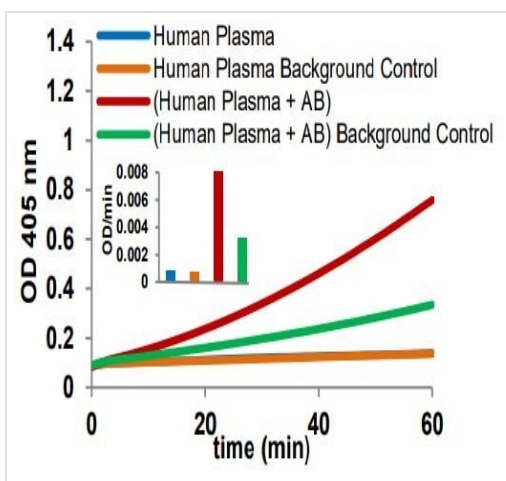


Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Measurement of purified Human Plasmin with or without the Plasmin Inhibitor Mix.

Example data



Measurement of Plasmin activity in Human Pooled EDTA-treated Plasma Sample (50 µl, 1:20 dilution) in the presence and absence of Activating Buffer containing Urokinase.

Example data

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