

Product datasheet

Urokinase type Plasminogen Activator Human Chromogenic Activity Assay Kit ab108915

[2 References](#) [3 Images](#)

Overview

Product name	Urokinase type Plasminogen Activator Human Chromogenic Activity Assay Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate
Assay type	Enzyme activity (quantitative)
Sensitivity	0.0051 U/ml
Species reactivity	Reacts with: Human
Product overview	<p>Urokinase type plasminogen activator (uPA) is a highly restricted serine protease that converts the zymogen plasminogen to active plasmin, a broad-spectrum serine proteinase capable of degrading most of the major protein components of the extracellular matrix. Binding of uPA to its receptor and subsequent uPA mediated pericellular proteolysis are involved in many process including cell migration and tissue remodelling in angiogenesis, atherogenesis, tumor cell metastasis, and ovulation. High level of uPA is a poor prognostic marker for aggressive breast cancer, aggressive prostate cancer, bladder cancer and gastric cancer.</p> <p>ab108915 uPA Human Chromogenic Activity Assay Kit is developed to determine Human uPA activity in plasma and cell culture supernatants. The assay measures the ability of uPA to activate the plasminogen to plasmin in coupled or indirect assays that contain uPA, plasminogen, and a plasmin-specific synthetic substrate. The amount of plasmin produced is quantitated using a highly specific plasmin substrate releasing a yellow para-nitroaniline (pNA) chromophore. The change in absorbance of the pNA in the reaction solution at 405 nm is directly proportional to the uPA enzymatic activity.</p> <p>The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.</p>
Platform	Microplate reader

Properties

Storage instructions

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

Components	100 tests
Assay Diluent	1 x 30ml
Human Plasminogen	1 x 2 vials
Human uPA Standard	1 vial
Microplate	1 unit
Plasmin Substrate	1 x 2 vials
Sealing Tapes	3 units

Function

Specifically cleave the zymogen plasminogen to form the active enzyme plasmin.

Tissue specificity

Expressed in the prostate gland and prostate cancers.

Sequence similarities

Belongs to the peptidase S1 family.

Contains 1 EGF-like domain.

Contains 1 kringle domain.

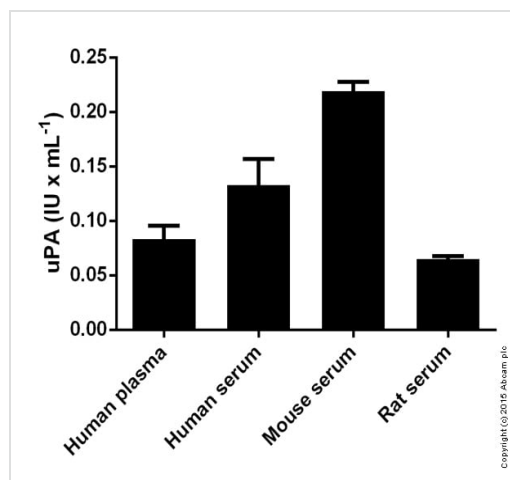
Contains 1 peptidase S1 domain.

Post-translational modifications

Phosphorylation of Ser-158 and Ser-323 abolishes proadhesive ability but does not interfere with receptor binding.

Cellular localization

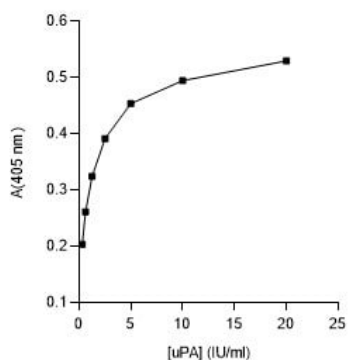
Secreted.

Images

uPA measured in biologicals showing quantity (IU) per mL of sample tested.

Indirect ELISA - Urokinase type plasminogen activator Human Chromogenic Activity Assay Kit (ab108915)

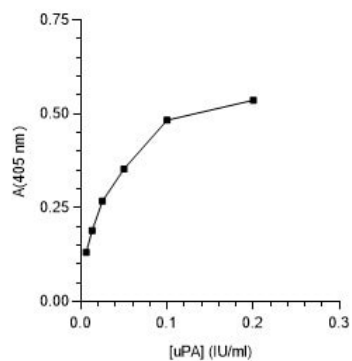
uPA Chromogenic Activity
Standard Curve (High uPA Activity Samples)



uPA Chromogenic Activity Standard Curve (For High uPA Activity Samples)

Indirect ELISA - Urokinase type plasminogen activator Human chromogenic activity assay kit
example standard curve (ab108915)

uPA Chromogenic Activity
Standard Curve (Low uPA Activity Samples)



uPA Chromogenic Activity Standard Curve (For Low uPA Activity Samples)

Indirect ELISA - Urokinase type plasminogen activator Human chromogenic activity assay kit
example standard curve (ab108915)

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