

# Human Urokinase type Plasminogen Activator ELISA Kit ab108917

[4 References](#) [1 Image](#)

### Overview

**Product name** Human Urokinase type Plasminogen Activator ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

| Sample  | n | Mean | SD | CV%  |
|---------|---|------|----|------|
| Overall |   |      |    | 4.9% |

Inter-assay

| Sample  | n | Mean | SD | CV%  |
|---------|---|------|----|------|
| Overall |   |      |    | 7.2% |

**Sample type** Cell culture supernatant, Urine, Serum, Plasma, Tissue

**Assay type** Sandwich (quantitative)

**Sensitivity** = 45 pg/ml

**Range** 0.125 ng/ml - 4 ng/ml

**Recovery** 98 %

**Assay time** 4h 00m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview**

Abcam's Urokinase type Plasminogen Activator Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Urokinase type Plasminogen Activator (uPA) in plasma, serum, urine, milk, cell culture supernatants and tissue extracts.

A Urokinase type Plasminogen Activator specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Urokinase type Plasminogen Activator specific biotinylated detection antibody is added and then followed by washing with wash buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize HRP enzymatic

reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Urokinase type Plasminogen Activator captured in plate.

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.**

**Platform** Microplate

## Properties

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

| Components   | 1 x 96 tests |
|--|--------------|
| 100X Streptavidin-Peroxidase Conjugate                               | 1 x 80µl     |
| 10X Diluent N Concentrate  | 1 x 30ml     |
| 20X Wash Buffer Concentrate  | 2 x 30ml     |
| 70X Biotinylated Human Urokinase type Plasminogen Activator Antibody | 1 x 120µl    |
| Chromogen Substrate  | 1 x 7ml      |
| Sealing Tapes  | 3 units      |
| Stop Solution  | 1 x 11ml     |
| Urokinase type Plasminogen Activator Microplate (12 x 8 well strips) | 1 unit       |
| Urokinase type Plasminogen Activator Standard                        | 1 vial       |

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

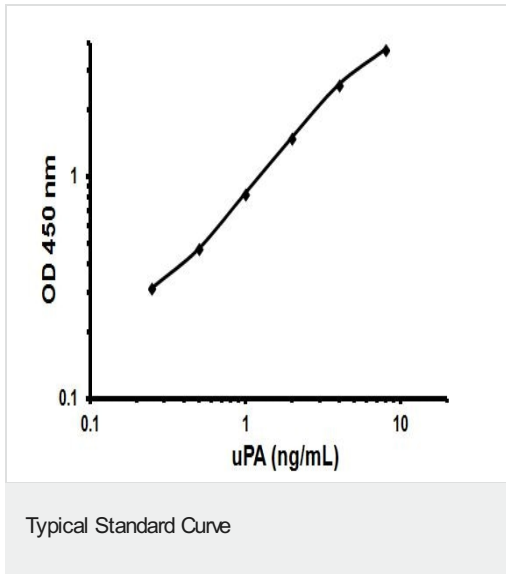
**Tissue specificity** Expressed in the prostate gland and prostate cancers.

**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



Representative Standard Curve using ab108917.

**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