

### Mouse uPA Total ELISA Kit (URK) ab198512

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

#### Overview

**Product name** Mouse uPA Total ELISA Kit (URK)

**Detection method** Colorimetric

#### Precision

Intra-assay

| Sample     | n  | Mean      | SD    | CV%   |
|------------|----|-----------|-------|-------|
| Known Conc | 20 | 0.06ng/ml | 0.005 | 8.12% |
| Known Conc | 20 | 0.3ng/ml  | 0.019 | 6.18% |
| Known Conc | 20 | 3.01ng/ml | 0.112 | 3.73% |

Inter-assay

| Sample     | n  | Mean      | SD    | CV%   |
|------------|----|-----------|-------|-------|
| Known Conc | 10 | 0.09ng/ml | 0.008 | 8.52% |
| Known Conc | 10 | 0.25ng/ml | 0.019 | 7.47% |
| Known Conc | 10 | 2.66ng/ml | 0.154 | 5.79% |

**Sample type** Plasma

**Assay type** Quantitative

**Sensitivity** 0.00998 ng/ml

**Range** 0.01 ng/ml - 10 ng/ml

**Recovery** 97.75 %

Sample specific recovery

| Sample type | Average % | Range        |
|-------------|-----------|--------------|
| Spike       | = 97.7    | 89.3% - 106% |

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Does not react with:** Sheep

## Product overview

Abcam's uPA Total (URK) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of total urokinase plasminogen activator antigen in mouse plasma.

Mouse uPA will bind to the capture antibody coated on the microtiter plate. Free and complexed enzyme will react with the capture antibody on the plate. After appropriate washing steps, polyclonal anti-murine uPA primary antibody binds to the captured enzyme. Excess antibody is washed away and bound polyclonal antibody is then bound by the secondary antibody conjugated to horseradish peroxidase. TMB substrate is used for color development at 450nm. A standard calibration curve is prepared along with the samples to be measured using dilutions of uPA.

## Platform

Microplate

## Properties

### Storage instructions

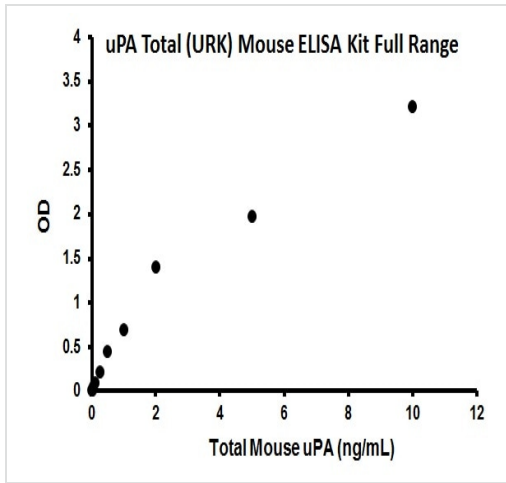
Store at +4°C. Please refer to protocols.

| Components  | 1 x 96 tests |
|---|--------------|
| 10X Wash Buffer   | 1 x 50ml     |
| Anti-mouse uPA primary antibody                                 | 1 vial       |
| Anti-rabbit horseradish peroxidase secondary antibody           | 1 vial       |
| Mouse uPA standard  | 1 vial       |
| TMB Substrate Solution  | 1 x 10ml     |
| uPA Antibody pre-coated 96 well microplate (12 x 8 well strips) | 1 x 96 tests |

## Relevance

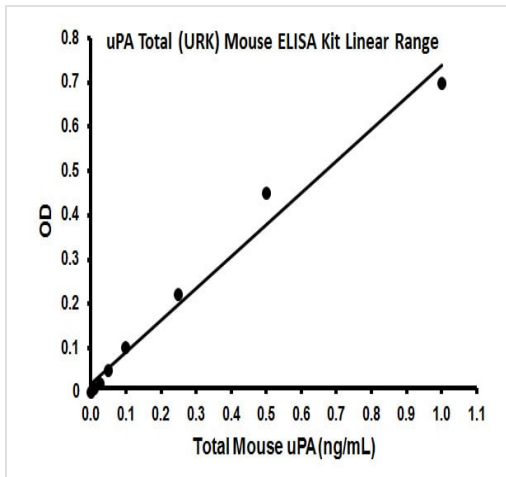
Urokinase type plasminogen activator (uPA) is secreted from cells as a precursor which is activated to the two chain form consisting of an A and B chain. The active high MW form is further processed by removal of an amino terminal fragment to an active low MW form (35 kDa). uPA is a serine protease that activates plasminogen to plasmin. High levels of uPA and plasminogen activator inhibitor type 1 (PAI 1) in breast cancer tissue extracts have been associated with rapid disease progression. The malignant phenotype of prostatic tumor cells correlates with the expression of both uPA and its cell membrane receptor (uPAR).

## Images



Full range standard curve using ab198512 uPA Total (URK) Mouse ELISA Kit.

Standard Curve.



Linear range standard curve using ab198512 uPA Total (URK) Mouse ELISA Kit.

Standard Curve.

**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