

Product datasheet

Mouse uPA ELISA Kit ab245727

Recombinant SimpleStep ELISA

[6 Images](#)

Overview

Product name Mouse uPA ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
serum	8			2.8%

Inter-assay

Sample	n	Mean	SD	CV%
serum	3			3.5%

Sample type Urine, Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 31.7 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Urine	112	101% - 124%
Serum	109	103% - 118%
Cell culture media	95	91% - 99%
Hep Plasma	110	105% - 114%
EDTA Plasma	113	108% - 120%
Cit plasma	106	104% - 106%

Assay time	1h 30m
Assay duration	One step assay
Species reactivity	Reacts with: Mouse Does not react with: Cow

Product overview

Mouse uPA ELISA Kit (ab245727) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of uPA protein in cit plasma, edta plasma, hep plasma, serum, urine, and cell culture media. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse uPA with 31.7 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Platform	Pre-coated microplate (12 x 8 well strips)
-----------------	--------------------------------------------

Properties

Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	-------------------------------------------

Components	1 x 96 tests
10X Mouse uPA Capture Antibody	1 x 600µl
10X Mouse uPA Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse uPA Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml

Components

1 x 96 tests

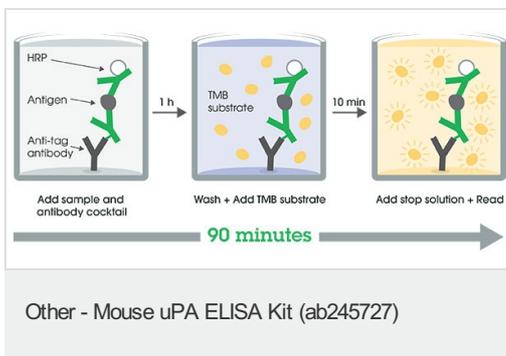
TMB Development Solution

1 x 12ml

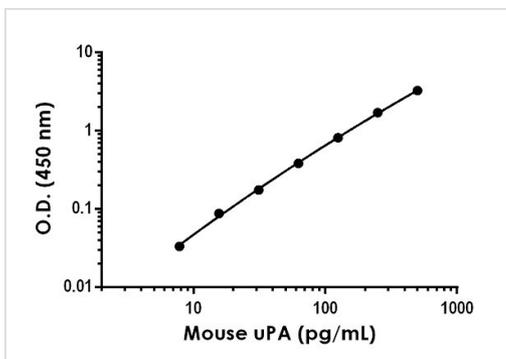
Relevance

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

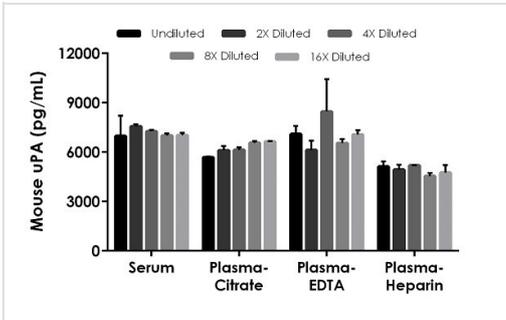


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



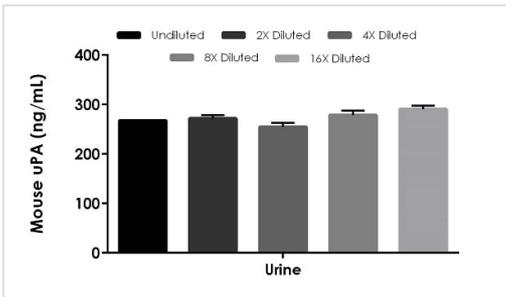
Example of mouse uPA standard curve in Sample Diluent NS.

The uPA standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



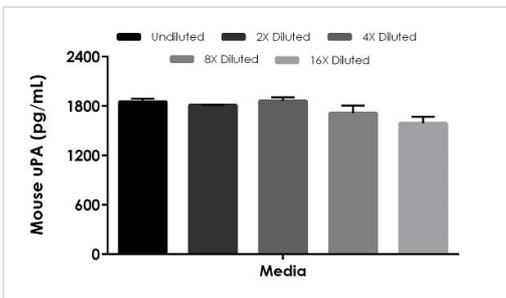
Interpolated concentrations of native uPA in mouse serum, plasma samples.

The concentrations of uPA were measured in duplicates, interpolated from the uPA standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (EDTA) 25%, and plasma (heparin) 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean uPA concentration was determined to be 7166 pg/mL in serum, 6222 pg/mL in plasma (citrate) and 7060 pg/mL in plasma (EDTA), and 4907 pg/mL in plasma (heparin).



Interpolated concentrations of native uPA in mouse urine.

The concentrations of uPA were measured in duplicates, interpolated from the uPA standard curves and corrected for sample dilution. Undiluted samples are as follows: urine 1.5%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean uPA concentration was determined to be 272.6 ng/mL in urine.



Interpolated concentrations of spiked uPA in cell culture media.

The concentrations of uPA were measured in duplicates, interpolated from the uPA standard curves and corrected for sample dilution. Undiluted samples are as follows: cell culture media 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Mouse uPA ELISA Kit
(ab245727)

To learn more about the advantages of recombinant antibodies see [here](#).

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