

Product datasheet

Uric Acid Assay Kit ab65344

[8 References](#) [4 Images](#)

Overview

Product name	Uric Acid Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Assay time	0h 40m
Product overview	Uric Acid Assay Kit (ab65344) provides a convenient means for detecting uric acid in biological samples such as serum and urine. Pretreatment of samples is not required.

Uric acid level can be measured using fluorometric (at Ex/Em = 535/587 nm) or colorimetric (at λ = 570 nm) methods.

Uric acid assay protocol summary:

- add reaction mix to samples and standards
- incubate for 30 min
- analyze on microplate reader

Notes Uric acid in serum is the end product of purine metabolism, and is cleared through the kidney by glomerular filtration. However, human often lacks the necessary enzyme called urate oxidase (Uricase), and therefore abnormal uric acid may be accumulated in blood. Recent evidences show the close association between serum urate level and cardiovascular morbidity and mortality, especially among persons at high cardiovascular risk, including those with hypertension, diabetes and congestion heart failure.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate reader

Properties

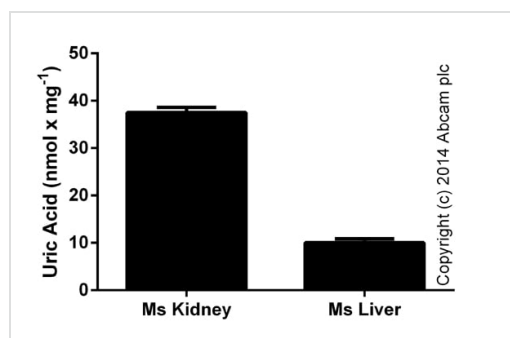
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	100 tests
Uric Acid Assay Buffer	WM	1 x 25ml
Uric Acid Enzyme Mix	Green	1 vial
Uric Acid Probe (in DMSO, anhydrous)	Red	1 x 0.2ml
Uric Acid Standard (2 nmol/μl)	Yellow	1 x 1ml

Relevance

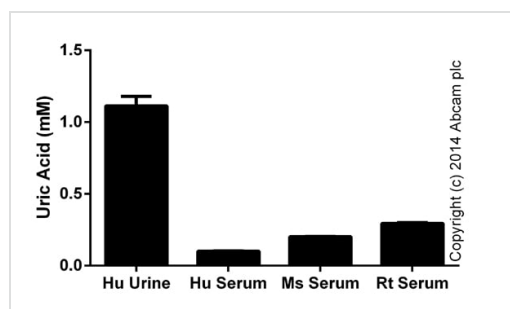
Uric acid is a heterocyclic purine derivate that is the final oxidation product of purine metabolism. It is produced by xanthine oxidase, which oxidizes oxypurines such as xanthine into uric acid. In most mammals, except humans and higher primates, the enzyme uricase further oxidizes uric acid to allantoin. In humans, however, uric acid is the end product of purine metabolism, and is cleared through the kidney by glomerular filtration and excreted in urine. Humans produce only small quantities of uric acid and therefore abnormal uric acid may be accumulated in blood leading to a type of arthritis known as gout.

Images



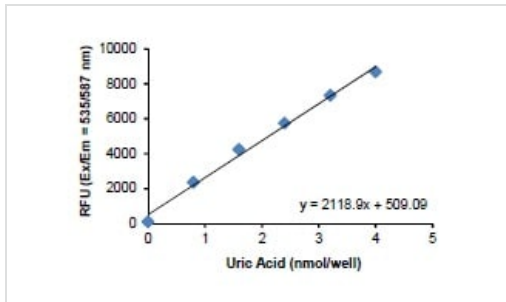
Uric acid measured in mouse tissue lysates, showing quantity (nmol) per mg of extracted protein (duplicates; +/- SD).

Functional assays: Uric acid (ab65344)



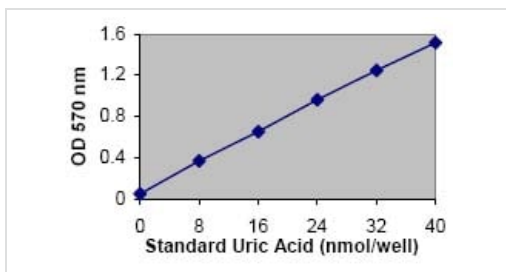
Uric acid measured in biological fluid (duplicates; +/- SD).

Functional assays: Uric acid (ab65344)



Uric acid fluorimetric standard curve.

Example of Fluometric Standard Curve



Uric acid colorimetric standard curve.

Functional Studies - Uric Acid Assay Kit (ab65344)

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