

Urea Assay Kit ab83362

★★★★★ 3 Abreviews 84 References 6 图像

概述

| | |
|------|--|
| 产品名称 | Urea Assay试剂盒 |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Urine, Serum, Plasma, Tissue Extracts, Cell Lysate |
| 检测类型 | Quantitative |
| 灵敏度 | > 0.5 nmol/well |
| 范围 | 0.5 nmol/well - 5 nmol/well |
| 检测时间 | 1h 00m |
| 产品概述 | <p>Urea Assay Kit ab83362 is a rapid, simple, sensitive, and reliable assay used to quantify urea in a variety of samples such as serum, plasma, and urine, etc. It can be used as a blood urea nitrogen assay kit (BUN assay kit).</p> <p>In the urea assay protocol, urea is acted on by enzymes to form a product that reacts with a probe to generate color (OD_{max}=570nm). The absorbance is directly proportional to the concentration of urea in the solution. The kit can detect as low as 0.5 nmol per well or 10 μM of urea.</p> <p>Urea assay protocol summary:</p> <ul style="list-style-type: none">- add samples and standards to wells- add reaction mix and incubate for 60 min at 37°C- analyze with microplate reader <p>This product is manufactured by BioVision, an Abcam company and was previously called K375 Urea Colorimetric Assay Kit. K375-100 is the same size as the 100 test size of ab83362.</p> <p>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.</p> <p>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.</p> |
| 平台 | Microplate reader |

性能

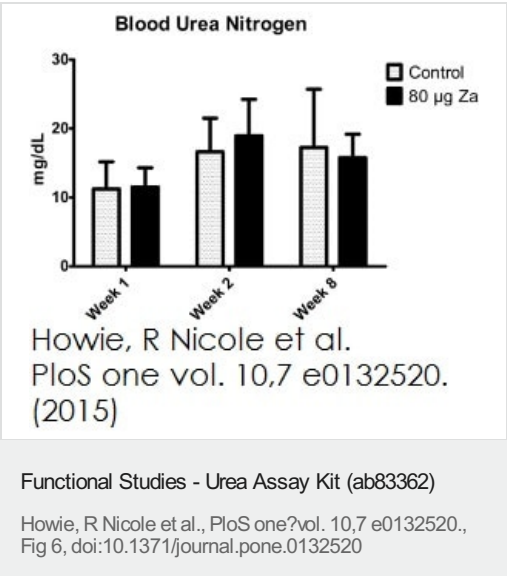
存放说明 Store at -20°C. Please refer to protocols.

| 组件 | 100 tests | 2000 tests |
|--------------------------|-----------|------------|
| Assay Buffer XIV | 1 x 25ml | 20 x 25ml |
| Converter Enzyme V | 1 vial | 20 vials |
| Developer I | 1 vial | 20 vials |
| Development Enzyme Mix I | 1 vial | 20 vials |
| OxiRed Probe | 1 x 200µl | 20 x 200µl |
| Urea Standard | 1 x 100µl | 20 x 100µl |

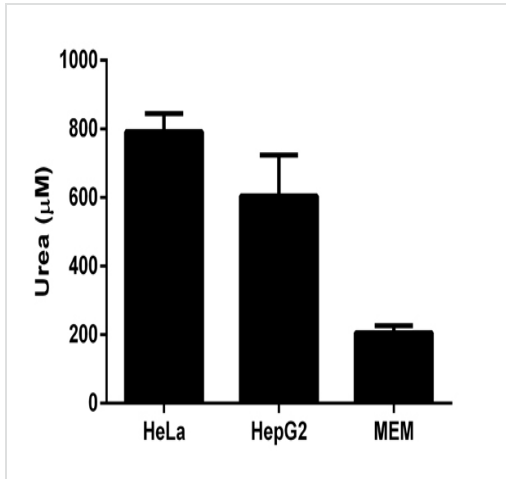
相关性

Urea is a waste product which is produced in the liver, dissolved in blood (in a concentration of 2.5 - 7.5 mM), and secreted by the kidneys. Urea also plays a very important role in protein catabolism, removal of toxic ammonia from the body, and the countercurrent system which allows for reabsorption of water and critical ions in the nephrons.

图片



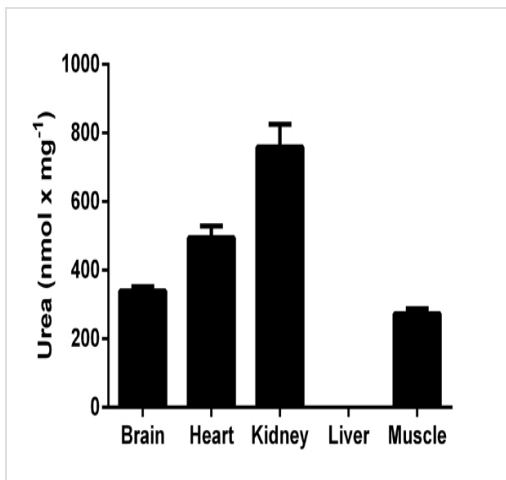
ab83362 was used as a BUN assay (Blood urea nitrogen assay) to measure urea levels in 80 female Sprague-Dawley rats (age 10-12 months-old).



Functional Studies - Urea Assay Kit (ab83362)

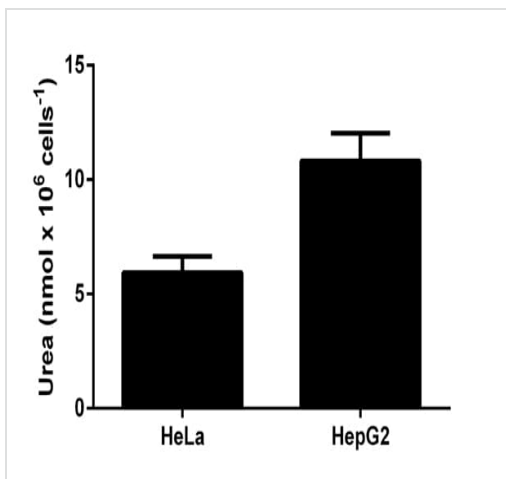
Urea measured in cell culture supernatants and control medium (MEM), background signal subtracted (duplicates +/- SD).

The samples were tested neat; 1:4; 1:16 and 1:64. Only the latter two gave reliable results.



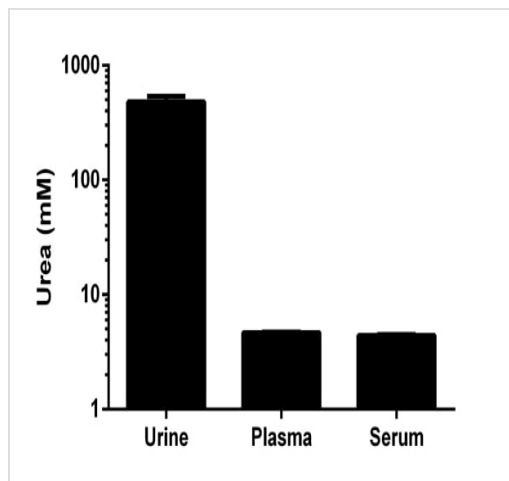
Functional Studies - Urea Assay Kit (ab83362)

Urea measured in mouse tissue lysates (mg of extracted protein), background signal subtracted (duplicates +/- SD).



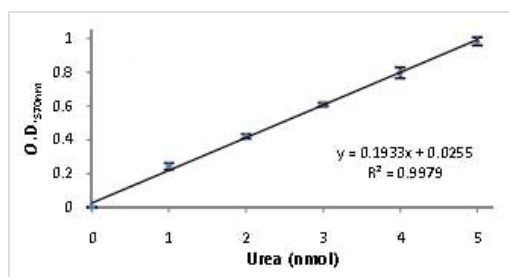
Functional Studies - Urea Assay Kit (ab83362)

Urea measured in cell lysates, background signal subtracted (duplicates; +/- SD).



Urea measured in biological fluids, background signal subtracted (duplicates; +/- SD).

Functional Studies - Urea Assay Kit (ab83362)



Example of Urea standard curve using ab83362

Functional Studies - Urea Assay Kit (ab83362)

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors