

Product datasheet

Creatinine Assay Kit ab65340

★★★★★ 2 Abreviews 25 References 5 Images

Overview

Product name	Creatinine Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids
Assay type	Quantitative
Assay time	1h 00m
Product overview	<p>Creatinine Assay Kit (ab65340) provides an accurate, convenient measure of creatinine concentration in biological fluids such as serum, urine or CSF. In the assay, creatinine is converted to creatine by creatininase, creatine is converted to sarcosine, which is specifically oxidized to produce a product which reacts with a probe to generate red color ($\lambda_{max} = 570 \text{ nm}$) and fluorescence ($Ex/Em = 538/587 \text{ nm}$). Unlike picric acid assays, this kit is suitable for serum/plasma creatinine determinations, as well as for urine and other biological samples.</p>

For deproteinization of samples: Better results are typically seen with this assay when using a 10kda filter for sample deproteinization than when using the PCA method.

Visit our [FAQs page](#) for tips and troubleshooting.

Notes	<p>Creatinine is a breakdown product of creatine phosphate. Creatinine is produced and excreted at a constant rate, and blood creatinine is used to determine glomerular filtration rate (GFR), a measure of kidney function. Blood creatinine levels increase only in cases of significant (>75%) damage to nephrons. Creatinine clearance is frequently used as a means of standardizing excretion of other compounds such as isoprostanes.</p>
--------------	--

Properties

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

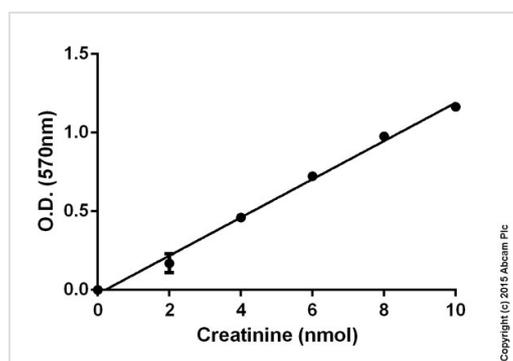
Components	Identifier	100 tests
Creatinase (Lyophilized)	Blue	1 vial
Creatininase (Lyophilized)	Violet	1 vial

Components	Identifier	100 tests
Creatinine Standard (10 μ mol) (Lyophilized)	Yellow	1 vial
Creatinine Assay Buffer	WM	1 x 25ml
Creatinine Enzyme Mix (Lyophilized)	Green	1 vial
Creatinine Probe in DMSO (200ul)		1 x 200 μ l

Relevance

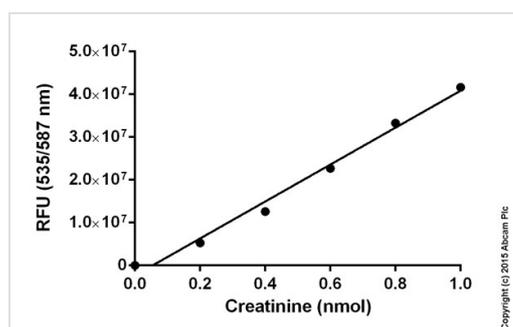
Creatinine, or creatine anhydride, is a breakdown product of creatine phosphate in muscle. The loss of water molecule from creatine results in the formation of creatinine. Creatinine is transferred to the kidneys by blood plasma, whereupon it is eliminated from the body by glomerular filtration and partial tubular excretion. Creatinine is usually produced and excreted at a fairly constant rate, and blood creatinine is used to determine glomerular filtration rate (GFR), a measure of kidney function.

Images



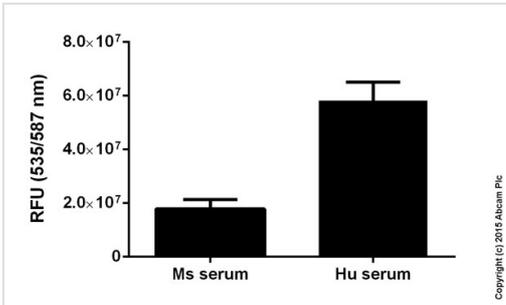
Standard curve: mean of duplicates (+/- SD)
with background reads subtracted

Functional Studies - Creatinine Assay Kit (ab65340)



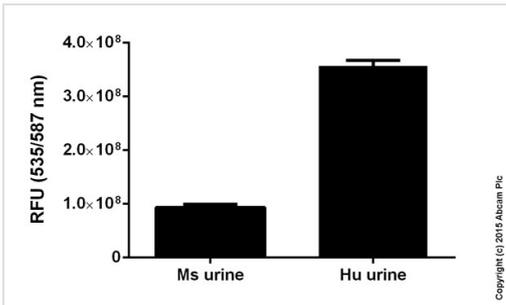
Standard curve: mean of duplicates (+/- SD)
with background reads subtracted

Functional Studies - Creatinine Assay Kit (ab65340)



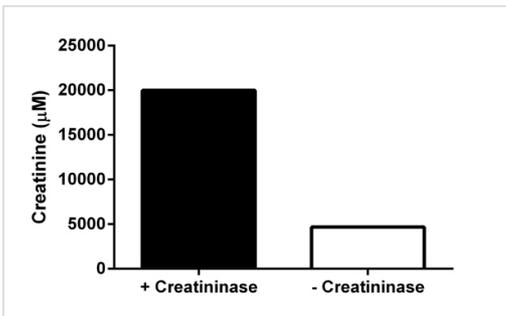
Creatinine measured in mouse and human serum plotted against RFU. Samples were diluted 5-10 fold.

Functional Studies - Creatinine Assay Kit (ab65340)



Creatinine measured in mouse and human urine plotted against RFU. Samples were diluted 400-800 fold.

Functional Studies - Creatinine Assay Kit (ab65340)



Creatinine levels in filtered human urine was measured in the presence or absence of creatinase (background signal subtracted).

Immunodiffusion - Creatinine Assay Kit (ab65340)

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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