

## Product datasheet

# Creatinine Assay Kit ab65340

★★★★★ 2 Abreviews 20 References 5 图像

### 概述

<b>产品名称</b>	Creatinine Assay试剂盒
<b>检测方法</b>	Colorimetric/Fluorometric
<b>样品类型</b>	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids
<b>检测类型</b>	Quantitative
<b>检测时间</b>	1h 00m
<b>产品概述</b>	<p>Creatinine Assay Kit (ab65340) provides an accurate, convenient measure of creatine 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 (<math>\lambda_{max} = 570 \text{ nm}</math>) and fluorescence (<math>E_x/E_m = 538/587 \text{ nm}</math>). Unlike picric acid assays, this kit is suitable for serum/plasma creatinine determinations, as well as for urine and other biological samples.</p> <p>For deproteinization of samples: 10kda method works the beter for deproteinization when preparing samples as opposed to PCA method. It's possible that the chemicals used negatively affects the enzyme components of the assay kit.</p>

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

<b>说明</b>	<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 (&gt;75%) damage to nephrons. Creatinine clearance is frequently used as a means of standardizing excretion of other compounds such as isoprostanes.</p>
-----------	--

<b>经测试应用</b>	<b>适用于:</b> Functional Studies
--------------	--------------------------------

### 性能

<b>存放说明</b>	Store at -20°C. Please refer to protocols.
-------------	--

组件	标识符	100 tests
Creatinase (Lyophilized)	Blue	1 vial

组件	标识符	100 tests
Creatininase (Lyophilized)	Violet	1 vial
Creatinine Standard (10 $\mu$ 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 $\mu$ l

**相关性** 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.

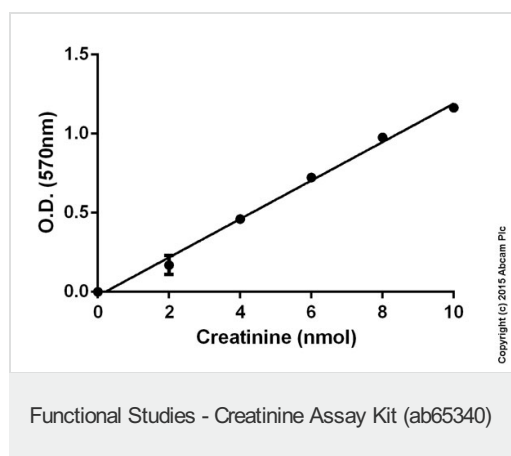
## 应用

Our [Abpromise guarantee](#) covers the use of **ab65340** in the following tested applications.

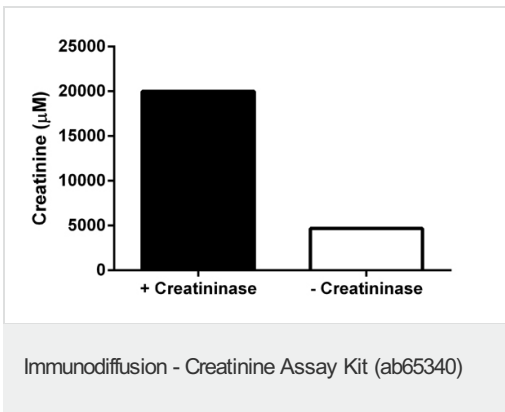
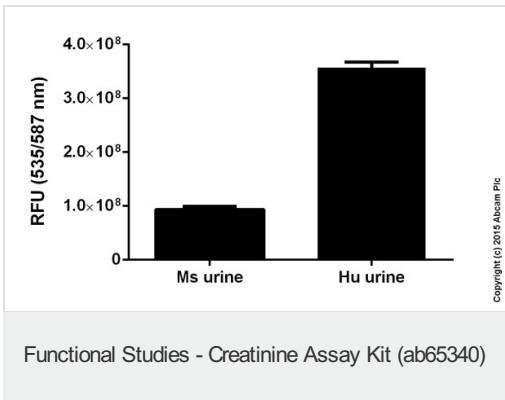
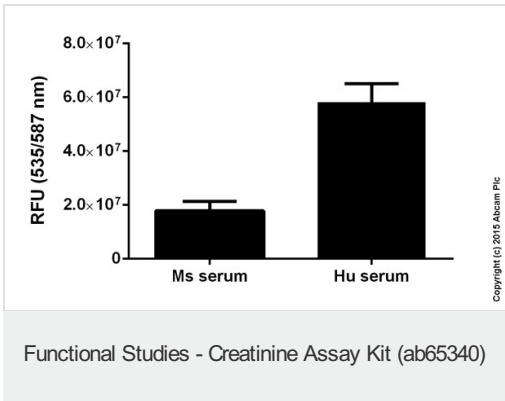
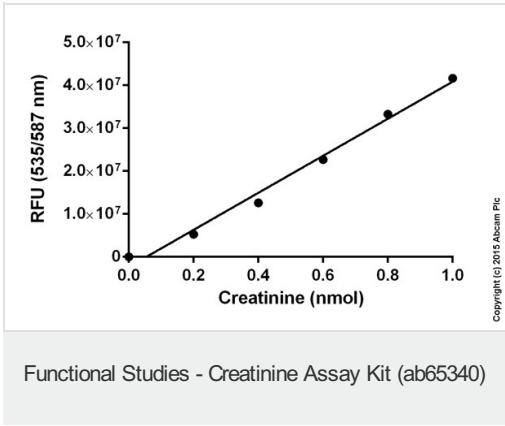
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Functional Studies		Use at an assay dependent dilution.

## 图片



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



**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 <http://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