

### Creatine Assay Kit (Colorimetric/Fluorometric) ab65339

[9 References](#) [2 图像](#)

#### 概述

产品名称	Creatine Assay试剂盒(Colorimetric/Fluorometric)
检测方法	Colorimetric/Fluorometric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Cell Lysate
检测类型	Quantitative
灵敏度	> 0.001 mM
范围	0.001 mM - 10 mM
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	<p>Creatine Assay Kit (Colorimetric/Fluorometric) (ab65339) provides an accurate, convenient measure of creatine in a variety of biological samples.</p> <p>In the creatine assay protocol, creatine is enzymatically converted to sarcosine which is then specifically oxidized to generate a product that converts a colorless probe to an intensely red color (<math>\lambda_{max} = 570 \text{ nm}</math>), and highly fluorescent (Ex/Em = 538/587 nm) product. Creatine is therefore easily detected by either colorimetric or fluorometric methods. Detection range 0.001–10 mM Creatine.</p> <p>Creatine assay protocol summary:</p> <ul style="list-style-type: none"><li>- add reaction mix to sample and standard wells</li><li>- incubate for 60 min</li><li>- analyze with a microplate reader</li></ul>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K635 Creatine Colorimetric/Fluorometric Assay Kit. K635-100 is the same size as the 100 test size of ab65339.</p> <p>Creatine is an endogenous compound whose function is to maintain a high ATP/ADP ratio, by way of its phosphorylated form and creatine kinase. Creatine supplementation has been used in the treatment of muscular, neurological and neurodegenerative diseases, as well as a sport performance enhancer. Detection of creatine level has wide applications in research and development.</p>
平台	Microplate reader

## 性能

## 存放说明

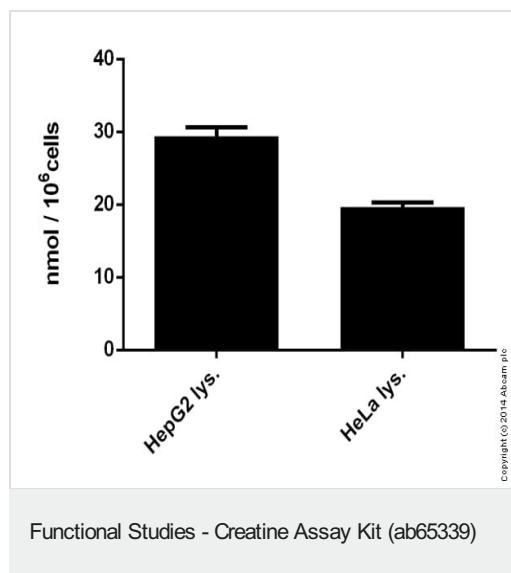
Store at -20°C. Please refer to protocols.

组件	100 tests
Assay Buffer II	1 x 25ml
Creatinase	1 vial
Creatine Standard	1 vial
OxiRed Probe	1 x 200µl
Sarcosine Enzyme Mix	1 vial

## 相关性

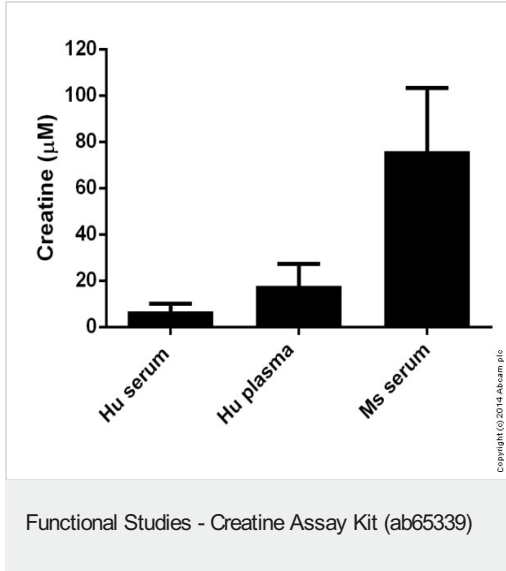
Creatine is an endogenous amino acid that occurs in vertebrate tissues and in urine. In muscle tissue, creatine is generally found as its phosphorylated form phosphocreatine. Creatine functions as part of the cell's energy shuttle: the high energy phosphate group of ATP is transferred to creatinine to form phosphocreatinine in the following reaction:  $Cr + ATP \rightarrow P-Cr + ADP$ . This reaction is reversibly catalyzed by creatine kinase. In the human body, creatine is synthesized mainly in the liver from arginine, glycine and methionine. Creatine is excreted as creatinine in the urine.

## 图片



Creatine measured in cell lysates showing quantity (nmol) per 10<sup>6</sup> cells

Functional Studies - Creatine Assay Kit (ab65339)



Creatine measured in biologicals showing concentration (micromolar)

**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