

### Creatinine Assay Kit (Colorimetric) ab204537

8 References [画像数 1](#)

#### 製品の概要

製品名 Creatinine Assay Kit (Colorimetric)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
7.981 mg/dl	20			= 2.159%
7.452 mg/dl	20			= 3.203%
2.007 mg/dl	20			= 5.995%
7.326 mg/dl	20			= 3.729%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
7.591 mg/dl	20			= 2.249%
7.232 mg/dl	20			= 2.052%
1.932 mg/dl	20			= 2.027%
7.085 mg/dl	20			= 1.433%

サンプルの種類 Urine

アッセイタイプ Quantitative

添加回収試験 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Urine	= 95	92% - 99%

全工程の試験時間 0h 30m

種交差性 交差種: Mammals, Other species

## 製品の概要

Abcam's Creatinine Assay Kit (Colorimetric) (ab204537) is a complete kit for the quantitative determination of creatinine in urine, and is based upon the Jaffe reaction.

**Sensitivity** = 0.042 mg/dL.

**General range** = 0.31 - 20 mg/dL

## 特記事項

Creatinine (2-amino-1-methyl-5H-imadazol-4-one) is a metabolite of phosphocreatine (p-creatine), a molecule used as a store for high-energy phosphate that can be utilized by tissues for the production of ATP. Creatine either comes from the diet or is synthesized from the amino acids arginine, glycine, and methionine. This occurs in the kidneys and liver, although other organ systems may be involved and species-specific differences may exist. Creatine and p-creatine are converted nonenzymatically to the metabolite creatinine, which diffuses into the blood and is excreted by the kidneys. In vivo, this conversion appears to be irreversible and in vitro it is favored by higher temperatures and lower pH2. Creatinine forms spontaneously from p-creatine, and under normal conditions, its formation occurs at a relatively constant rate. Intra-individual variation of creatinine levels is <15% from day to day, making it a useful marker for normalizing levels of other molecules found in urine. Altered creatinine levels may be associated with conditions that result in decreased renal blood flow, such as diabetes and cardiovascular disease.

## 試験プラットフォーム

Microplate reader

## 製品の特性

### 保存方法

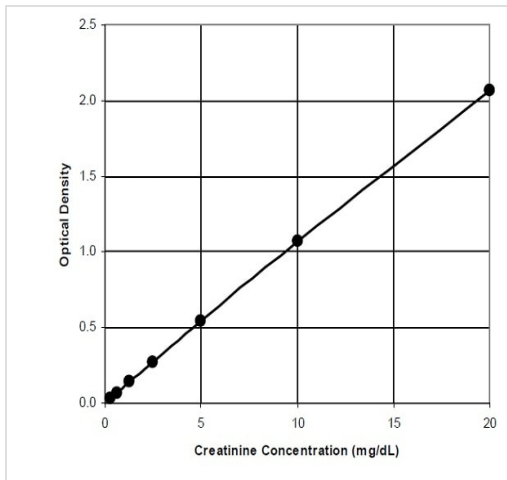
Store at +4°C. Please refer to protocols.

内容	2 x 96 tests
Clear Microtiter Plates	2 units
Creatinine Detection Reagent	1 x 20ml
Creatinine Standard	1 vial
Plate Sealer	2 units

## 関連性

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.

## 画像



Typical standard curve obtained using Creatinine Assay Kit (Colorimetric) (ab204537) which is a complete kit for the quantitative determination of creatinine in urine, and is based upon the Jaffe reaction.

Typical standard curve obtained using Creatinine Assay Kit (Colorimetric) (ab204537).

**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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

---

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