

Product datasheet

Protein Creatinine Ratio Assay Kit ab272539

2 Images

Overview

Product name	Protein Creatinine Ratio Assay Kit
Detection method	Colorimetric
Sample type	Urine
Assay type	Quantitative
Product overview	<p>Protein Creatinine Ratio Assay Kit (ab272539) is a simple and convenient test for proteinuria. Other methods such as 24 hour urine test or timed urine test require strict adherence to sample collection protocol. Since the protein concentration is normalized to creatinine secretion, the urine sample can be taken at anytime and no diet or liquid restrictions are necessary for sample collection.</p> <p>Sensitive and accurate: Use 20 µL samples. Linear detection range in 96-well plate: 1 - 20 mg/dL Protein and 1 – 150 mg/dL Creatinine.</p> <p>Fast and convenient: No sample pre-treatment is needed. Simple 10- minute "add-incubate-read" procedure.</p> <p>High-throughput adaptable: The procedure can be readily automated for processing thousands of samples per day.</p>
Tested applications	Suitable for: Functional Studies

Properties

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

Components	100 tests
CR Reagent A	1 x 6ml
CR Reagent B	1 x 6ml
PR Reagent	1 x 24ml
Standard	1 x 1ml

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.

Applications

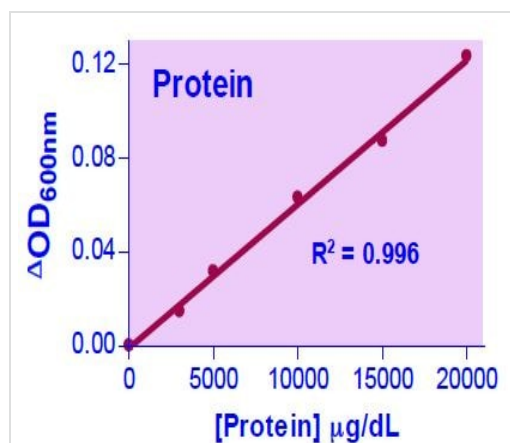
The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab272539 in the following tested applications.

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

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

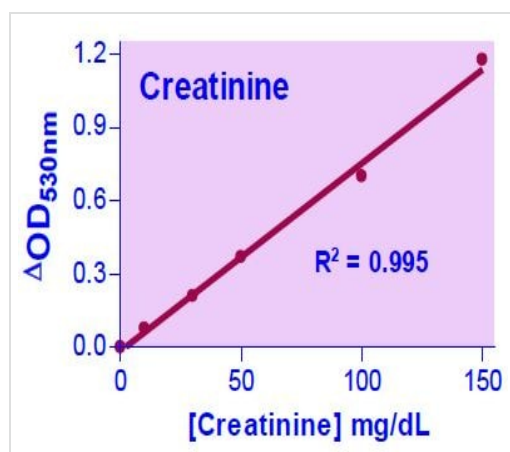
Images



Protein Standard Curve

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Example Protein standard curve.



Creatinine Standard Curve

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Example Creatinine standard curve.

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.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