

Coenzyme A Assay Kit ab102504

[6 References](#) [2 图像](#)

概述

产品名称	Coenzyme A Assay试剂盒
样品类型	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
检测类型	Quantitative
灵敏度	> 2.5 μ M
范围	2.5 μ M - 250 μ M
检测时间	1h 00m
产品概述	Abcam's Coenzyme A Assay Kit provides an easy, accurate assay to measure the CoA level in variety biological samples. In the assay, free CoA is specifically utilized to generate products which react with OxiRed Probe to generate color ($\lambda = 570$ nm) and fluorescence (Ex=535/Em=587 nm). The assay can detect 0.1 to 10 nmol of CoA (2.5-250 μ M concentration range) in a variety of samples.

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

说明 This product is manufactured by BioVision, an Abcam company and was previously called K367 Coenzyme A (CoA) Colorimetric/Fluorometric Assay Kit. K367-100 is the same size as the 100 test size of ab102504.

Coenzyme A (CoA) is composed of units derived from cysteine, pantothenic acid, and ATP. It plays important roles in the synthesis and oxidation of fatty acids, pyruvate oxidation in the citric acid cycle and many other biological processes. One of the main functions of CoA is the carrying and transfer of acyl groups. One of the most important acyl groups transferred is the acetate group, in which case the molecule is called acetyl-CoA. The acetyl group eventually finds itself incorporated into a variety of molecules such as cholesterol, acetylcholine, melatonin, heme and the TCA cycle intermediates.

平台 Microplate reader

性能

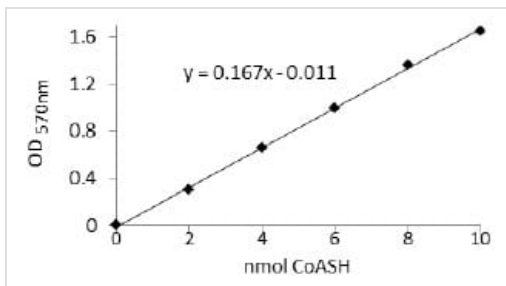
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
ACS Substrate	1 x 1ml

组件	100 tests
Acyl CoA Enzyme Mix	1 vial
Assay Buffer V	1 x 25ml
CoA Converter Mix	1 unit
CoA Standard	1 unit
OxiRed Probe	1 unit

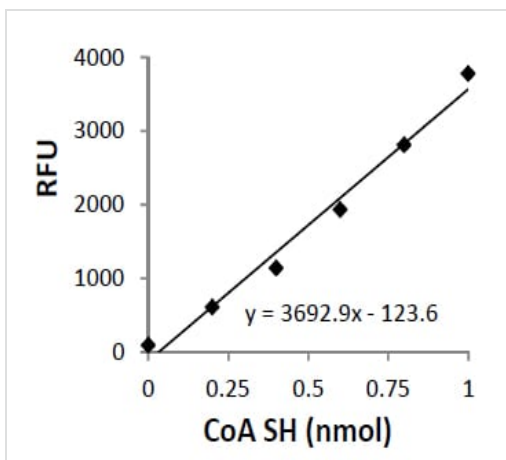
相关性 Coenzyme A (CoA) is composed of units derived from cysteine, pantothenic acid, and ATP. It plays important roles in the synthesis and oxidation of fatty acids, pyruvate oxidation in the citric acid cycle and many other biological processes. One of the main functions of Coenzyme A is the carrying and transfer of acyl groups. One of the most important acyl groups transferred is the acetate group, in which case the molecule is called acetyl-Coenzyme A. The acetyl group eventually finds itself incorporated into a variety of molecules such as cholesterol, acetylcholine, melatonin, heme and the TCA cycle intermediates.

图片



Example of Standard Curve obtained using ab102504.

Functional Studies - Coenzyme A Detection Kit (ab102504)



Example of Standard Curve obtained using ab102504.

Functional Studies - Coenzyme A Detection Kit (ab102504)

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