

### Coenzyme A Assay Kit (Fluorometric - Green) ab138889

6 References 画像数 1

#### 製品の概要

製品名	Coenzyme A Assay Kit (Fluorometric - Green)
検出方法	Fluorescent
サンプルの種類	Cell culture extracts, Tissue Extracts, Cell Lysate
アッセイタイプ	Quantitative
検出感度	= 40 nM
全工程の試験時間	1h 00m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Coenzyme A Assay Kit (Fluorometric - Green) (ab138889) provides an ultrasensitive fluorometric assay to quantitate CoA content by detection of the –SH group in CoA.</p> <p>The fluorogenic CoA green indicator dye used in the kit becomes strongly fluorescent upon reacting with –SH.</p> <p>The assay kit can detect as little as 4 pmol CoA/ 100 µL assay volume (40 nM). It can be performed in a convenient 96-well or 384-well microtiter-plate format at Ex/Em = 490/520 nm, and easily adapted to automation without a separation step.</p> <p>Coenzyme A assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add samples and standards to wells</li> <li>- add reaction mix</li> <li>- incubate for 10-60 min whilst measuring fluorescence with a microplate reader</li> </ul> <p><b>The kit is for solution-based inhibitors or activators of Coenzyme A only. Any samples that contain proteins or thiol-containing compounds are not compatible with the kit.</b></p>
特記事項	Previously called Coenzyme A Detection kit (Fluorometric - Green).
試験プラットフォーム	Microplate reader

#### 製品の特性

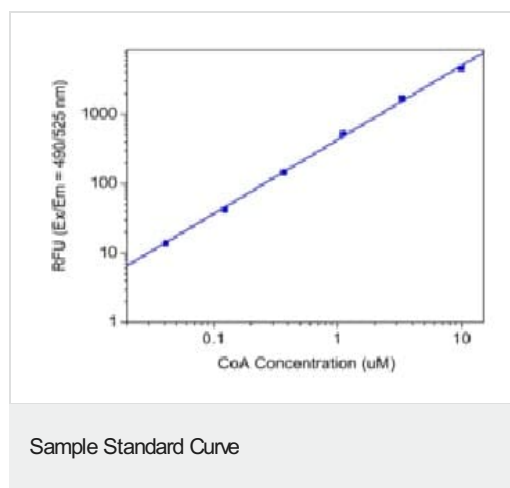
保存方法 Store at -20°C. Please refer to protocols.

内容	200 tests
CoA Green Indicator	1 vial
Assay Buffer	1 x 25ml
Coenzyme A Standard	1 x 154µg
DMSO	1 x 200µl

**関連性**

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.

## 画像



CoA dose response was measured in a 96-well black plate with Coenzyme A Detection Kit using a microplate reader. As low as 40 nM (4 pmol/well) of CoA was detected with 30 minutes incubation time (n=3).

**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