

Isocitrate Assay Kit ab83424

画像数 1

製品の概要

製品名	Isocitrate Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell Lysate, Tissue Lysate, Food samples
アッセイタイプ	Quantitative
検出感度	> 0.2 µg
検出範囲	0.2 µg - 5 µg
全工程の試験時間	0h 40m
製品の概要	Isocitrate Assay Kit (ab83424) provides a simple, sensitive and rapid means of quantifying isocitrate in a variety of samples. In this assay, isocitrate is oxidized with the generation of NADPH which converts a nearly colorless probe to an intensely colored species with a λmax of 450nm. This assay can detect 1 to 20 nmoles (~0.2-5 µg) of isocitrate.

	Visit our <a href="#">FAQs page</a> for tips and troubleshooting.
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K656 Isocitrate Colorimetric Assay Kit. K656-100 is the same size as the 100 test size of ab83424.</p> <p>Isocitric acid (HOOC-CHOH-CH (-COOH)-CH2-COOH) is an intermediate of the Krebs TCA cycle, positioned between citrate and α-keto-glutarate. It is the branch point from which the glyoxylate shunt operates in plants and lower organisms. Isocitrate is found in substantial concentrations in many fruits and vegetables as well as in foods produced from these raw materials. In the TCA cycle, isocitrate is oxidized by isocitrate dehydrogenase (IDH) to α-ketoglutarate with the generation of NAD(P)H. Loss of NAD-IDH has been implicated as a potential causative factor in retinitis pigmentosa.</p>
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.
------	--

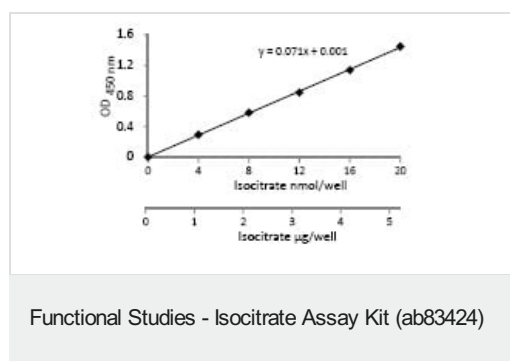
内容	100 tests
Developer (Lyophilised)	1 vial

内容	100 tests
Isocitrate Assay Buffer	1 x 25ml
Isocitrate Enzyme Mix	1 x 200µl
Isocitrate Standard	1 x 100µl

## 関連性

Isocitric acid (HOOC-CHOH-CH (-COOH)-CH<sub>2</sub>-COOH) is an intermediate of the Krebs TCA cycle, positioned between citrate and alpha-keto-glutarate. It is the branch point from which the glyoxylate shunt operates in plants and lower organisms. Isocitrate is found in substantial concentrations in many fruits and vegetables as well as in foods produced from these raw materials. In the TCA cycle, isocitrate is oxidized by isocitrate dehydrogenase (IDH) to alpha-ketoglutarate with the generation of NAD(P)H. Loss of NAD-IDH has been implicated as a potential causative factor in retinitis pigmentosa.

## 画像



Isocitrate standard curve generated using this kit protocol.

**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