

Malate Assay Kit ab83391

5 References 画像数 1

製品の概要

製品名	Malate Assay Kit
検出方法	Colorimetric
サンプルの種類	Tissue Extracts, Food samples
アッセイタイプ	Quantitative
検出感度	20 µM
検出範囲	20 µM - 200 µM
全工程の試験時間	1h 00m
製品の概要	Abcam's Malate Assay Kit is an easy and sensitive assay to measure the L(-) Malate level in a variety of samples. In the assay, malate is specifically oxidized to generate a product which reacts with a substrate probe to generate color (λmax = 450 nm). The assay can detect 1~10 nmol of Malate in a 50 µl sample with a detection sensitivity ~20 µM.

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

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K637 Malate Colorimetric Assay Kit. K637-100 is the same size as the 100 test size of ab83391.</p> <p>L(-) Malate is a TCA cycle intermediate. It plays an important role in the Calvin cycle during carbon fixation in plants. In lower organisms, malate is converted to lactate during malolactic fermentation with the formation of CO2. Malate is frequently used as an additive in the food and pharmaceutical industries, so quantitating malic acid is important in manufacturing beer, wine, cheese and fruits, among others.</p>
------	---

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

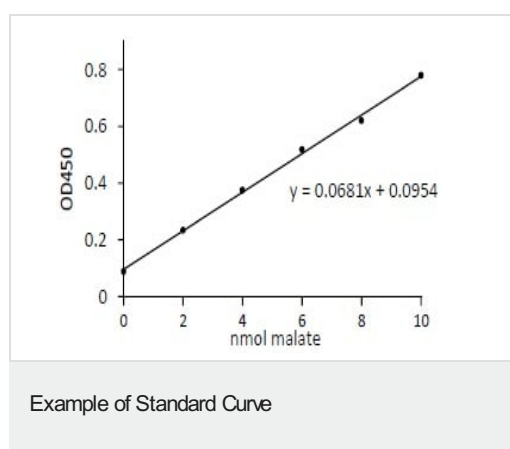
保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	100 tests	100 tests
Assay Buffer LXIV	1 x 25ml	1 x 25ml
Developer Solution III	1 vial	1 vial

内容	100 tests	100 tests
Malate Enzyme Mix	1 vial	1 vial
Malate Standard	1 vial	1 vial

関連性 L(-) Malate is a TCA cycle intermediate. It plays an important role in the Calvin cycle during carbon fixation in plants. In lower organisms, malate is converted to lactate during malolactic fermentation with the formation of CO₂. Malate is frequently used as an additive in the food and pharmaceutical industries, so quantitating malic acid is important in manufacturing beer, wine, cheese and fruits, among others.

画像



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