

Malate Dehydrogenase Activity Assay Kit (Colorimetric) ab183305

3 References **画像数 3**

製品の概要

製品名	Malate Dehydrogenase Activity Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Enzyme activity (quantitative)
検出感度	< 0.5 mU/well
種交差性	交差種: Mammals, Other species
製品の概要	In Abcam's Malate Dehydrogenase Activity Assay kit (Colorimetric) (ab183305), malate dehydrogenase (MDH) reacts with malate to form an intermediate, which in turn reacts with MDH Developer to form a colored product with strong absorbance at 450 nm. The assay is simple, sensitive and can detect less than 0.5 mU of MDH activity in various sample types.

This assay detects both isoforms of MDH: MDH1 and MDH2. If you want to detect enzymatic activity of each specific isoform, please consider **MDH1 Activity Assay (ab200009)** or **MDH2 activity Assay (ab119693)**.

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K654 Malate Dehydrogenase Activity Colorimetric Assay Kit. K654-100 is the same size as the 100 test size of ab183305.

Malate Dehydrogenase (MDH) (EC 1.1.1.37) is an important enzyme which reversibly converts L-malate into oxaloacetate in the presence of NAD. In eukaryotic cells, malate dehydrogenase has 2 isoforms: MDH1 and MDH2. MDH1 is cytosolic & participates in the malateaspartate shuttle, which transports malate into mitochondria for utilization in ATP generation whereas MDH2 is a mitochondrial enzyme and part of the citric acid cycle. MDH activity is increased in some neurodegenerative diseases such as Alzheimer's disease, and abnormal MDH activity in serum can serve as a diagnostic tool for severe liver damage (e.g. Hepatocellular carcinoma).

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Microplate reader

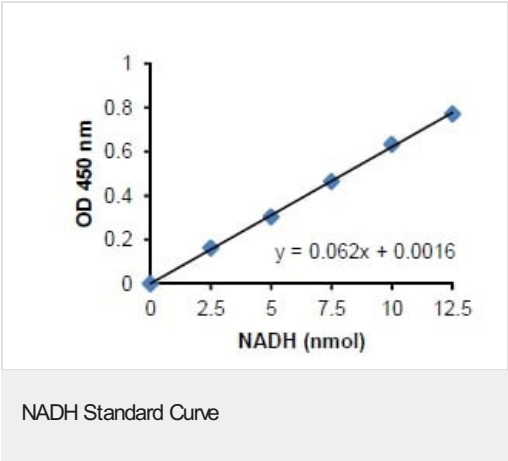
製品の特性

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

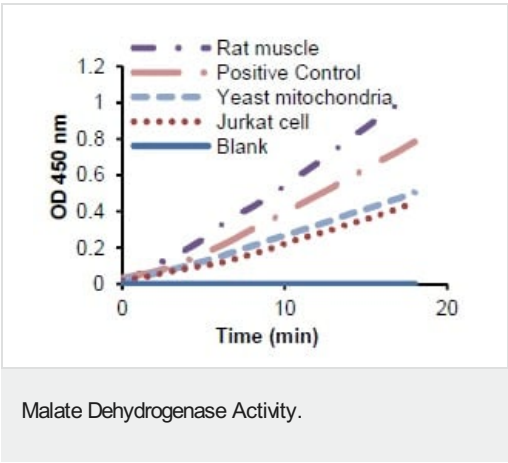
内容	100 tests
MDH Assay Buffer	1 x 25ml
MDH Developer (Lyophilized)	1 vial
MDH Enzyme Mix (Lyophilized)	1 vial
MDH Positive Control (Lyophilized)	1 vial
MDH Substrate (Lyophilized)	1 vial
NADH Standard (Lyophilized)	1 vial

細胞内局在MDH2: Mitochondrion matrix. MDH1: Cytoplasm.

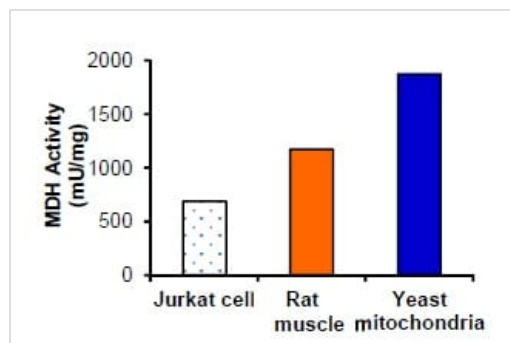
画像



NADH Standard Curve. Malate Dehydrogenase activity in rat muscle extract (0.8 µg), Jurkat cell lysate (0.6 µg), yeast mitochondria (1.2 µg) & MDH Positive Control.



Rat muscle extract (0.8 µg)
Jurkat cell lysate (0.6 µg)
Yeast mitochondria (1.2 µg)
MDH positive control.



Referenced MDH Activity in Jurkat cell lysate, rat muscle extract and yeast mitochondrial lysate.

Referenced MDH activity in Jurkat cell lysate, rat muscle extract and yeast mitochondrial lysate.

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