

α-Ketoglutarate Dehydrogenase Activity Assay Kit (Colorimetric) ab185440

[13 References](#) [3 Images](#)

Overview

Product name	a-Ketoglutarate Dehydrogenase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Semi-quantitative
Sensitivity	< 100 μU
Product overview	Abcam's α-Ketoglutarate Dehydrogenase Activity Assay kit (Colorimetric) (ab185440) provides a quick and easy way for monitoring α-KGDH activity in various samples. In the assay, α-KGDH converts α-ketoglutarate into an intermediate which reduces the probe to a colored product with strong absorbance at 450 nm. The assay is simple, sensitive and can detect α-ketoglutarate dehydrogenase activity lower than 0.1 mU in a variety of samples.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K678 Alpha-Ketoglutarate Dehydrogenase Activity Colorimetric Assay Kit. K678-100 is the same size as the 100 test size of ab185440.</p> <p>α-Ketoglutarate Dehydrogenase (α-KGDH) (EC 1.2.4.2) is a key enzyme in the citric acid cycle. It forms an enzyme complex with dihydrolipoamide succinyl transferase (E2) and dihydrolipoamide dehydrogenase (E3). α-KGDH converts α-ketoglutarate into succinylCoA in the presence of NAD and CoA. It is highly regulated by intracellular ATP/ADP and NADH/NAD ratios and calcium. In humans, decreased KGDH activity can lead to neurodegenerative diseases such as Alzheimer's disease. Recent studies show that α-KGDH is a target of oxidative stress; reactive oxygen species (ROS) inhibit KGDH activity which diminishes its critical function and can cause a bioenergetic deficit.</p>
Platform	Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer IX	1 x 25ml

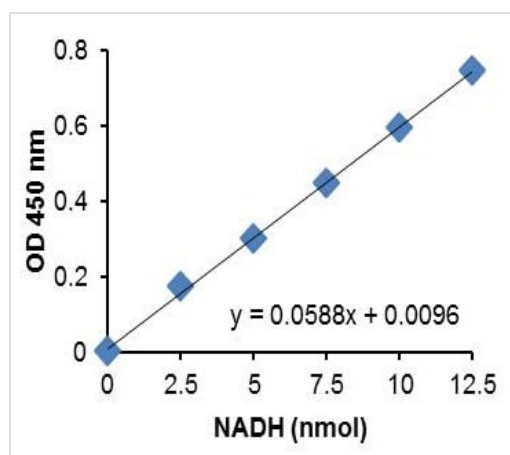
Components	100 tests
Developer Solution III	1 vial
KGDH Positive Control	1 x 50µl
KGDH Substrate Mix	1 vial
NADH Standard I	1 vial

Function The 2-oxoglutarate dehydrogenase complex catalyzes the overall conversion of 2-oxoglutarate to succinyl-CoA and CO₂. It contains multiple copies of three enzymatic components: 2-oxoglutarate dehydrogenase (E1), dihydrolipoamide succinyltransferase (E2) and lipoamide dehydrogenase (E3).

Sequence similarities Belongs to the alpha-ketoglutarate dehydrogenase family.

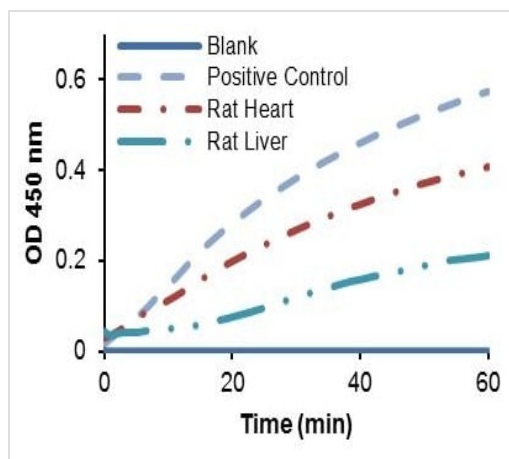
Cellular localization Mitochondrion matrix.

Images



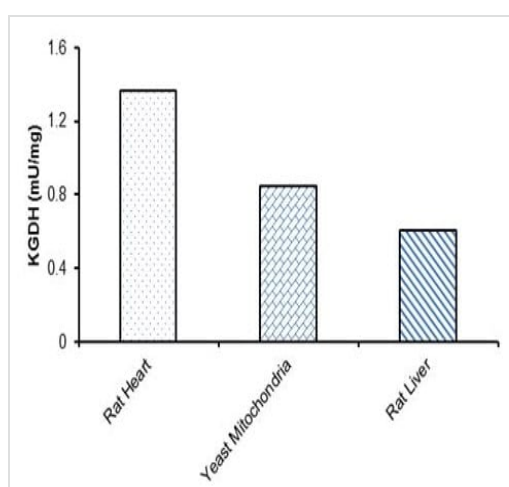
This is example data only.

NADH Standard Curve



alpha-Ketoglutarate Dehydrogenase activity in various samples

α -Ketoglutarate Dehydrogenase activity in rat heart (75 μ g) and liver lysates (100 μ g). Assays were performed following the kit protocol. This is example data only.



alpha-Ketoglutarate Dehydrogenase specific activity in various samples

α -Ketoglutarate Dehydrogenase specific activity was calculated in rat heart lysate (75 μ g), yeast mitochondria prepared from *S. Cerevisiae* (10 μ g) and in rat liver lysate (100 μ g). Assays were performed following the kit protocol. This is example data only.

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors