

Succinate Assay Kit (Colorimetric) ab204718

★★★★★ [1 Abreviews](#) [36 References](#) [画像数 2](#)

医薬用外毒物

製品の概要

製品名	Succinate 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	Quantitative
検出感度	40 µM
全工程の試験時間	0h 30m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Succinate Assay Kit (Colorimetric) (ab204718) is a sensitive, fast and easy-to-use kit. In the succinate assay protocol, Succinate is utilized by Succinyl-CoA Synthetase to form an intermediate, which undergoes a series of reactions and reduces a colorless probe to a colored product with strong absorbance at OD=450 nm. This assay kit can detect less than 40 µM of succinate or succinic acid in a variety of sample types.</p> <p>Succinate + ATP + CoA converted by succinyl-CoA synthase to succinyl-CoA + ADP + Pi with Color Detection (OD 450 nm).</p> <p>Succinate assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix and incubate for 30 min - analyze with a microplate reader
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K649 Succinate (Succinic Acid) Colorimetric Assay Kit. K649-100 is the same size as the 100 test size of ab204718.</p> <p>Succinic acid (C₄H₆O₄) is distributed in all plants and animal tissues and was first obtained from amber. Due to its low toxicity, it is widely used in agriculture, food and pharmaceutical industry. Succinate (C₄H₄O₄) is the salt or ester of succinic acid. It is one of the most active components of cellular respiratory and intracellular energy generation. Measurement of succinate or succinic acid level is a key to analysis of the citric acid cycle.</p> <p>This kit has so far only been tested with human and rat samples. Since Succinate is distributed in all plants and animal tissues this kit should detect Succinate independent of the species. Depending on the sample type, the sample preparation might need to be optimized.</p>

製品の特性

保存方法

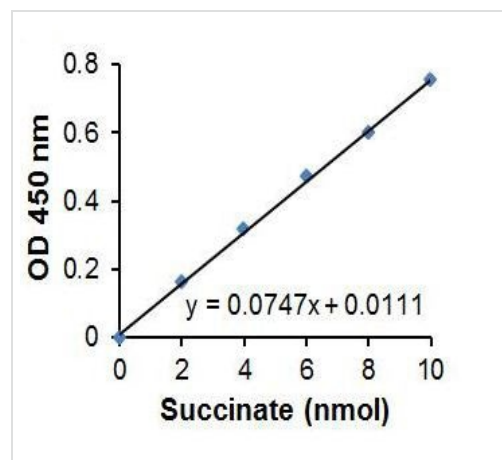
Store at -20°C. Please refer to protocols.

内容	100 tests
Developer Solution III	1 vial
Enzyme Mix XI	1 vial
Assay Buffer VII	1 x 25ml
Succinate Converter (Lyophilized)	1 vial
Succinate Standard	1 vial
Succinate Substrate Mix	1 vial

関連性

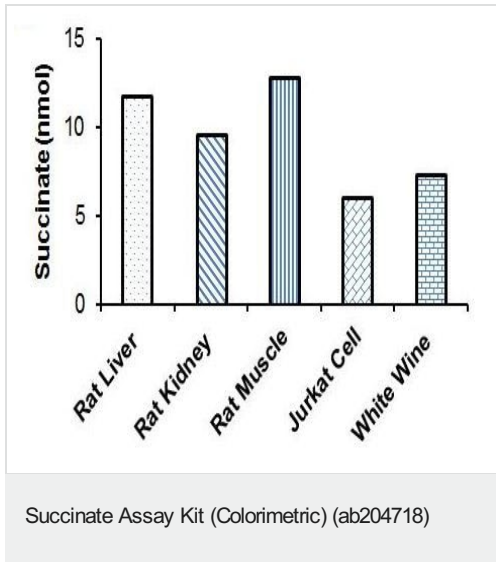
Succinic acid is involved in the catabolism of the neurotransmitter gamma-aminobutyric acid (GABA).

画像



Succinate Standard Curve.

Succinate Assay Kit (Colorimetric) (ab204718)



Measurement of Succinate levels in rat liver (40 µg), kidney (40 µg) and muscle (40 µg), Jurkat cell lysate (200 µg) and white wine (4 µl, wine sample was diluted 1:1 with 0.5 M Tris HCl, pH 8.0).

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