

Citrate Synthase Activity Assay Kit ab119692

★★★★★ [1 Abreviews](#) [28 References](#) [画像数 2](#)

製品の概要

製品名	Citrate Synthase Activity Assay Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	1	3		CV%
				5.45%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	1	3		CV%
				8.3%
サンプルの種類	Cell culture extracts, Tissue Extracts			
アッセイタイプ	Enzyme activity			
検出感度	1.56 µg/ml			
検出範囲	1.56 µg/ml - 100 µg/ml			
全工程の試験時間	3h 15m			
種交差性	交差種: Cow, Human 非交差種: Mouse, Rat			
製品の概要	<p>Citrate Synthase Activity Assay Kit (ab119692) is used to determine mitochondrial citrate synthase (CS) activity in a sample. The enzyme is immunocaptured within the wells of a microplate, and activity is determined by recording color development of TNB, which is generated in the reaction of citrate synthesis. The reaction product, TNB, absorbs at OD = 412 nm.</p> <p>This kit immunocaptures only native citrate synthase from the samples. Using the immunocapture method allows the measurement of citrate synthase activity with simple sample preparation and no need for mitochondrial isolation.</p> <p>This kit can be used with whole cell extracts or tissue homogenates.</p> <p>Citrate synthase assay protocol summary:</p> <ul style="list-style-type: none">- add samples to wells			

- incubate for 3 hrs to capture enzyme
- wash with wash buffer
- add activity solution
- analyze with microplate reader for 5-30 min at 20s intervals

特記事項

Citrate synthase (CS, O75390) is the initial enzyme of the tricarboxylic acid (TCA) cycle. This enzyme is a 51.7 kDa enzyme that catalyzes the reaction of 2 carbon acetyl CoA with 4 carbon oxaloacetate to form the 6 carbon citrate (EC 2.3.3.1). This enzyme is an exclusive marker of the mitochondrial matrix. Citrate synthase is found in nearly all cells capable of oxidative metabolism.

Storage: All components are shipped cold. 10X Reaction Reagent mixture is shipped lyophilized. Before use, rehydrate the reaction mixture by adding 1mL base buffer. Vortex the tube thoroughly to dissolve the lyophilized powder to prepare 10X Reaction solution. After hydration unused amounts of the 10X Reaction solution should be stored at -80°C for 6 months. Store all other components at 4°C. This kit is stable for 6 months from receipt.

試験プラットフォーム

Microplate reader

製品の特性

保存方法

Please refer to protocols.

内容	1 x 96 tests
10X Blocking Buffer	1 x 6ml
10X Reaction Reagent	1 vial
1X Base Buffer	1 x 25ml
20X Buffer	1 x 20ml
CS Microplate	1 unit
Extraction Buffer (ab260490)	1 x 15ml

パスウェイ

Carbohydrate metabolism; tricarboxylic acid cycle; isocitrate from oxaloacetate: step 1/2.

配列類似性

Belongs to the citrate synthase family.

細胞内局在

Mitochondrion matrix.

画像

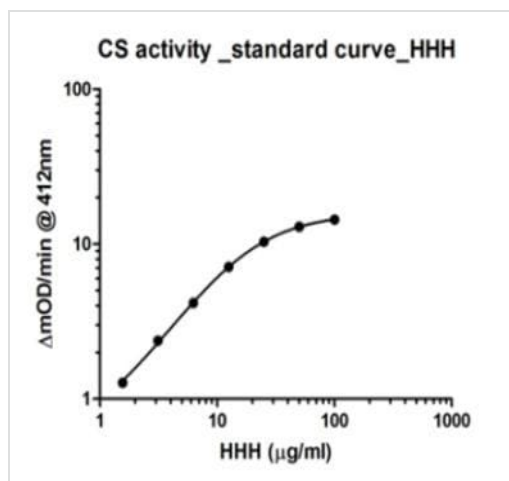


Figure 2. Example activity measurements from titrated of Human heart homogenate.

ELISA - Citrate Synthase Activity Assay Kit
(ab119692)

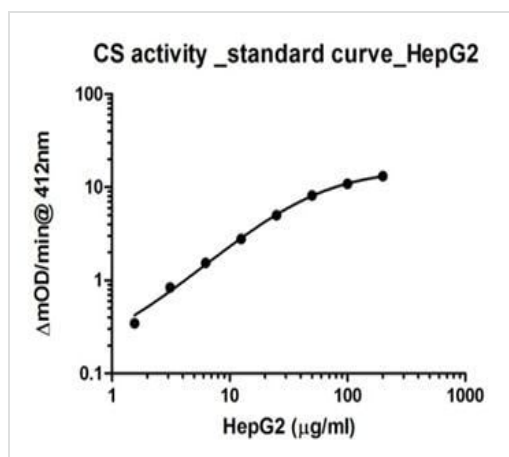


Figure 1. Example activity measurements from titrated HepG2 cells.

ELISA - Citrate Synthase Activity Assay Kit
(ab119692)

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