



ab119692

Human Citrate Synthase (CS) Activity Assay Kit

Instructions for Use

For the measurement of mitochondrial citrate synthase (CS) activity in human

This product is for research use only and is not intended for diagnostic use.

www.abcam.com

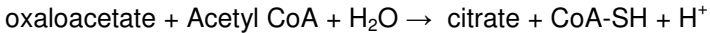
ab119692 Citrate Synthase Activity Assay Kit

Table of Contents

1. Introduction	3
2. Assay Summary	6
3. Kit Contents	7
4. Storage and Handling	8
5. Additional Materials Required	8
6. Reagent Preparation	9
7. Test Sample Preparation	12
8. Assay Procedure	14
9. Data Analysis	19

1. Introduction

Principle: The microplate assay ab119692 is used to determine mitochondrial citrate synthase (CS) activity in a sample at an immunocapture based manner. The enzyme is captured within the wells of the microplate, and activity is determined by recording color development of TNB, which is generated from DTNB present in the reaction of citrate synthesis. The overall reaction product, TNB, absorbs at 412 nm.



(↑ Absorbance at 412 nm)

ab119692 specifically immunocaptures only native citrate synthase in well from the applied test sample. The immunocapture based activity assay allows researchers to measure citrate synthase activity with simple sample preparation and no need for mitochondria isolation.

The kit contains all the required reagents for a fast and simple measurement of citrate synthase activity in a whole cell extract or in tissue homogenate.

Background: 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.

Limitations:

- FOR RESEARCH US ONLY. NOT FOR DIAGNOSTIC PROCEDURES.
- Use this kit before expiration date.
- Do not mix or substitute reagents from other lots or sources.
- If samples generate values outside of the range of the standard curve, further dilute the samples with 1X Incubation buffer and repeat the assay.
- Any variation in operator, pipetting technique, washing technique, incubation time or temperature, and kit age can cause variation in binding.

Technical Hints:

- To avoid cross contamination, change pipette tips between each addition of standard, sample and other reagents. Also use separate clean, dry reservoirs for each reagent.
- Cover plate during incubation steps.
- Thorough and consistent wash technique is essential for proper assay performance. Wash buffer must be forcefully dispensed and completely removed from the wells by aspiration or decanting. Remove remaining wash buffer by inverting the plate and blotting on paper towels.

2. Assay Summary

Prepare samples as instructed.

Determine the protein concentration of extracts.



Bring all reagents to room temperature.



Dilute sample to desired protein concentration in incubation buffer.

Add 100 μ L sample to each well used. Incubate 3 hours at room temperature.



Aspirate and wash each well twice.



Freshly prepare 1X Activity Solution. Add 100 μ L 1X Activity Solution to each well.

Pop bubbles and record immediately the color development with time at 412 nm for 15 to 30 minutes.

3. Kit Contents

Sufficient materials are provided for 96 measurements in a microplate.

Item	Quantity
20X Buffer	20 ml
Extraction Buffer	15 ml
10X Blocking Buffer	6 ml
Base Buffer	25 mL
10X Reaction Reagent	1 Vial
CS Microplate (12 x 8 antibody coated well strips)	96 Wells

4. Storage and Handling

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.

5. Additional Materials Required

- Standard absorbance microplate reader capable of kinetic reading at 412nm
- Multichannel pipette (50 - 300 μ L) and tips
- 1.5-mL microtubes
- Paper towels
- Deionized water
- Optional plate shaker for all incubation steps

6. Reagent Preparation

- 6.1. Prepare 1X Wash Buffer by adding 20 mL 20X Buffer to 380 mL nanopure water.
- 6.2. Prepare 1X Incubation Buffer by adding 6 mL 10X Blocking Solution to 54 mL 1X Wash Buffer.
- 6.3. Prepare 10X Reaction Solution, as described above (4. storage and handling).
- 6.4. Before use (in Step 8.6) prepare 1X Activity Solution. For the entire plate add 9mL Base Buffer to 1mL 10X Reaction Solution.

Note – the final concentration of substrate is now 1mM oxaloacetate, 0.06mM Acetyl CoA, and 0.1mM DTNB.

7. Test sample Preparation

Note: Extraction buffer DOES NOT CONTAIN PROTEASE INHIBITORS. It can be supplemented with phosphatase inhibitors, PMSF and protease inhibitor cocktail prior to use.

Supplements should be used according to manufacturer's instructions.

7.1 Cell lysates or extracts

- 7.1.1. Collect non adherent cells by centrifugation or scrape to collect adherent cells from the culture flask. Typical centrifugation conditions for cells are 500 x g for 10 min at 4°C.
- 7.1.2. Rinse cells twice with PBS.
- 7.1.3. Solubilize cell pellet at 2×10^7 /mL in Extraction Buffer.
- 7.1.4. Incubate on ice for 20 minutes. Centrifuge at 16000 x g 4°C for 20 minutes. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C for 6 months. The sample protein concentration in the extract may be quantified using a protein assay.

7.2 Tissue lysates

- 7.2.1. Tissue lysates are typically prepared by homogenization of tissue that is first minced and

ab119692 Citrate Synthase Activity Assay Kit

thoroughly rinsed in PBS to remove blood (dounce homogenizer recommended).

- 7.2.2. Suspend the homogenate to 25 mg/mL in PBS.
- 7.2.3. Solubilize the homogenate by adding 4 volumes of Extraction Buffer to a sample protein concentration of 5 mg/mL.
- 7.2.4. Incubate on ice for 20 minutes. Centrifuge at 16,000 x g for 20 minutes at 4 °C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80 °C for 6 months. The sample protein concentration in the extract may be quantified using a protein assay.

7.3 Sub-cellular organelle lysates e.g. mitochondria:

- 7.3.1. Prepare the organelle sample by, for example, sub-cellular fractionation.
- 7.3.2. Pellet the sample.
- 7.3.3. Solubilize the pellet by adding 9 volumes Extraction Buffer.

ab119692 Citrate Synthase Activity Assay Kit

- 7.3.4. Incubate on ice for 20 minutes. Centrifuge at 16,000 x g for 20 minutes at 4 °C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80 °C for 6 months. The sample protein concentration in the extract may be quantified using a protein assay.

These test samples should be diluted to within the working range of the assay in 1X Incubation Buffer.

8. Assay Procedure

Equilibrate all reagents and samples to room temperature before use. It is recommended all samples and standards be assayed in duplicate.

- 8.1. Prepare all reagents, and samples as directed in the previous sections.
- 8.2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, and re-seal.
- 8.3. Add 100 μ L of each diluted sample per well. It is recommended to include a dilution series of a control

ab119692 Citrate Synthase Activity Assay Kit

(normal) sample as a reference. Also include a 1X Incubation Buffer as a zero standard.

- 8.4. Cover/seal the plate and incubate for 3 hours at room temperature. If available use a plate shaker for all incubation steps at 300 rpm.
- 8.5. Aspirate each well and wash, repeat this once more for a total of two washes. Wash by aspirating or decanting from wells then dispensing 300 μ L 1X Wash buffer into each well as described above. Complete removal of liquid at each step is essential to good performance. After the last wash, remove the remaining buffer by aspiration or decanting. Invert the plate and blot it against clean paper towels to remove excess liquid.
- 8.6. Gently add 100 μ L fresh mixed 1X Activity Solution, described above, to each well minimizing the production of bubbles.
- 8.7. Pop any bubbles immediately and record absorbance in the microplate reader prepared as follows:

Mode:	Kinetic
Wavelength:	412 nM
Time:	5 - 30 min.
Temperature	Ambient
Interval:	20 sec.
Shaking:	Shake between readings

- 8.8. *Assay reactions occur faster as ambient temperatures are increased.
- 8.9. Analyze the data as described below.

9. Data Analysis

Example data sets are shown below illustrating data analysis of CS activity measurements in HepG2 cells (as an example human cell line derived from liver), and homogenate samples from human, rat and mouse liver tissues.

The starting concentration of Oxaloacetate and Acetyl CoA in the assay is 1mM and 0.06mM respectively. DTNB is added to the citrate synthase reaction to produce TNB, which can be measured at 412nm. For simplicity the activity can be expressed as the change in absorbance per minute per amount of sample loaded into the well. Activity was collected as described in this protocol using a Molecular Dynamics microplate reader. Standard curves of reference sample data were exported to graphing software capable of a 4-parameter

data analysis (shown below). Activity is clearly measurable in the 1.56-100 $\mu\text{g}/\text{mL}$ range when such a fit is applied for HepG2, and human tissue extracts.

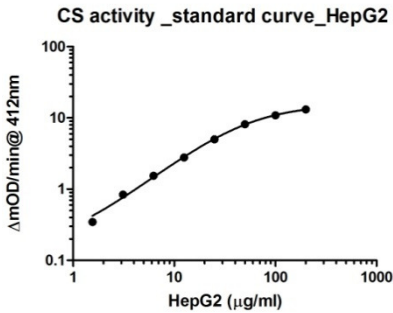


Figure 1. Example activity measurements from titrated HepG2 cells.

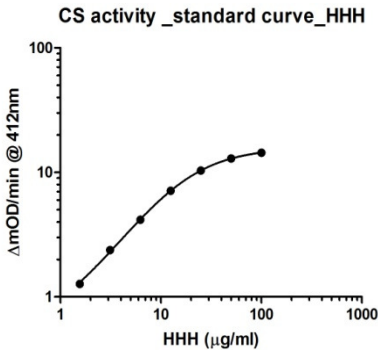


Figure 2. Example activity measurements from titrated of Human heart homogenate.

Unknown samples should be interpolated from these reference sample graphs. This determined relative activity is the amount of reference sample required to generate the same amount of activity as the unknown sample and usually expressed as a percent value.

WORKING RANGE

This assay has been demonstrated with human, rat, and mouse liver and heart tissue homogenate samples as well as HepG2 whole cell lysate, a liver derived human cultured cell line. Typical ranges for several sample types are described below. It is highly recommended to prepare multiple dilutions for each sample to ensure that each is in the working range of the assay (see Data Analysis section).

Sample Type	Range %
Cultured whole cell extracts (type dependant) e.g. HepG2	1.56 - 100 µg/mL
Tissue extract e.g. heart homogenate	1.56 - 100 µg/mL

REPRODUCIBILITY

Parameter	%CV
Intra (n=3, 3 assays at 2,525 µg/mL)	4.35-6.55
Inter (n=3 days, at 25 µg/mL))	8.3

SPECIFICITY

1. Species– human. Does not react with rat and mouse. Others not tested.

ab119692 Citrate Synthase Activity Assay Kit

2. The antibody used to isolate CS in this kit was generated by immunization of bovine heart mitochondrial matrix. The resulting monoclonal mouse antibody specifically immunoprecipitates a single CS band to purity from human samples. The immunoprecipitate was confirmed to be CS by mass spectrometry with no other contamination. The specificity of the immunocapture based activity assay is confirmed by replacing the capture antibody with another unrelated capture antibody.

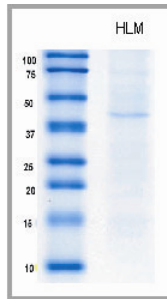


Figure 3. Monoclonal mouse antibody from this kit immunoprecipitated with human liver mitochondria exhibiting a single CS band.

CS activity _specificity_ HepG2/H3@25 μ g/mL

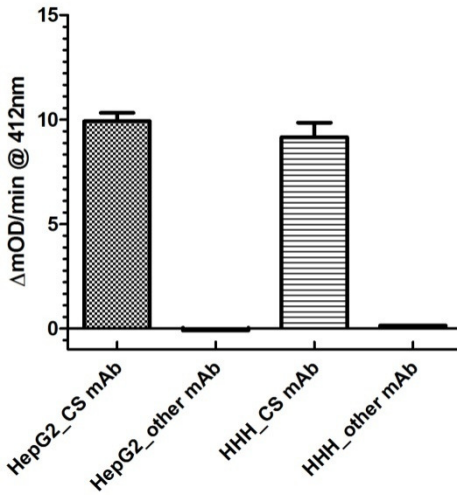


Figure 4. Immunocapture based activity assay specificity is confirmed by replacing the capture antibody with another unrelated capture antibody.

11. Troubleshooting

Problem	Cause	Solution
Poor standard curve	Inaccurate Pipetting	Check pipets
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Low cleaved PARP concentration in sample	Use appropriate positive control/treatment to induce apoptosis/PARP cleavage
	Incubation times too brief	Ensure sufficient incubation times; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions
	Contaminated wash buffer	Make fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Store your reconstituted standards at -80 °C, all other assay components 4 °C. Keep substrate solution protected from light

ab119692 Citrate Synthase Activity Assay Kit

ab119692 Citrate Synthase Activity Assay Kit

ab119692 Citrate Synthase Activity Assay Kit

Abcam in the USA

Abcam Inc
1 Kendall Square, Ste B2304
Cambridge,
MA 02139-1517
USA

Toll free: 888-77-ABCAM (22226)
Fax: 866-739-9884

Abcam in Europe

Abcam plc
330 Cambridge Science Park
Cambridge
CB4 0FL
UK

Tel: +44 (0)1223 696000
Fax: +44 (0)1223 771600

Abcam in Japan

Abcam KK
2-2-1 Nihonbashi
Horidome-cho,
Chuo-ku, Tokyo
103-0012
Japan

Tel: +81-(0)3-6231-094
Fax: +81-(0)3-6231-0941

Abcam in Hong Kong

Abcam (Hong Kong) Ltd
Unit 225A & 225B, 2/F
Core Building 2
1 Science Park West Avenue
Hong Kong Science Park
Hong Kong

Tel: (852) 2603-682
Fax: (852) 3016-1888