

Product Information

Succinyl-CoA Synthetase Activity Colorimetric Assay Kit

Catalog Number **MAK217**
Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

Succinyl-CoA synthetase (SCS; succinyl-CoA ligase; succinate thiokinase; EC 6.2.1.5) is a mitochondrial enzyme that hydrolyzes ATP to convert succinate to succinyl-CoA in the tricarboxylic acid (Krebs) cycle.¹ Deficiency in SCS activity is associated with encephalomyopathy and depletion of mitochondrial DNA.²

The Succinyl-CoA Synthetase Activity Colorimetric Assay Kit is a simple and high throughput adaptable assay for measuring SCS activity in a variety of tissues, cells, and isolated mitochondria. SCS activity is determined by measuring a colorimetric product with absorbance at 450 nm (A_{450}) proportional to the enzymatic activity present. One unit of succinyl-CoA synthetase is the amount of enzyme that generates 1.0 μmole of NADH per minute at pH 7.4 at $25\text{ }^{\circ}\text{C}$.

Components

The kit is sufficient for 100 assays in 96 well plates.

| | |
|--|-------|
| SCS Assay Buffer Catalog Number MAK217A | 25 mL |
| SCS Substrate Mix Catalog Number MAK217B | 1 vL |
| SCS Enzyme Mix Catalog Number MAK217C | 1 vL |
| SCS Developer Catalog Number MAK217D | 1 vL |
| NADH Standard Catalog Number MAK217E | 1 vL |
| SCS Positive Control Catalog Number MAK217F | 1 vL |

Reagents and Equipment Required but Not Provided

- 96 well flat-bottom plate – It is recommended to use clear plates for colorimetric assays.
- Spectrophotometric multiwell plate reader
- Mitochondria Isolation Kit (optional for mitochondria samples, Catalog Number MITOISO1 for tissue, MITOISO2 for cells, MITOISO3 for yeast, or equivalent)

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Briefly centrifuge vials before opening. Use ultrapure water for the preparation of reagents. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

SCS Assay Buffer – Store the buffer at $-20\text{ }^{\circ}\text{C}$ or $2-8\text{ }^{\circ}\text{C}$. Allow buffer to come to room temperature before use.

SCS Substrate Mix and SCS Developer – Reconstitute with 220 μL of water. Dissolve completely and mix by pipetting. Aliquot and store at $-20\text{ }^{\circ}\text{C}$. Keep on ice during use. Use within 2 months.

SDH Enzyme Mix – Reconstitute with 220 μL of SCS Assay Buffer. Dissolve completely and mix by pipetting. Aliquot and store at $-20\text{ }^{\circ}\text{C}$. Keep on ice during use. Use within 2 months.

NADH Standard – Reconstitute with 400 μL of water to generate a 1.25 mM NADH Standard Solution. Aliquot and store at $-20\text{ }^{\circ}\text{C}$. Keep on ice during use. Use within 2 months.

SCS Positive Control – Reconstitute with 100 μL of SCS Assay Buffer. Mix by pipetting. Aliquot and store at $-70\text{ }^{\circ}\text{C}$. Keep on ice during use. Use within 2 months.

Storage/Stability

The kit is shipped on wet ice and storage at $-20\text{ }^{\circ}\text{C}$, protected from light, is recommended.

Procedure

All samples and standards should be run in duplicate.

NADH Standards

Add 0, 2, 4, 6, 8, and 10 μL of the 1.25 mM (1.25 nmole/ μL) NADH Standard Solution into a 96 well plate, generating 0 (blank), 2.5, 5.0, 7.5, 10, and 12.5 nmole/well standards. Add SCS Assay Buffer to each well to bring the volume to 50 μL .

Sample Preparation

Homogenize tissue (10 mg) or cells (1×10^6) in 100 μL of ice cold SCS Assay Buffer. Keep on ice for 10 minutes. Centrifuge at $10,000 \times g$ for 5 minutes at $4\text{ }^{\circ}\text{C}$ minutes to remove insoluble material. Transfer supernatant to a fresh tube.

When analyzing SCS activity in mitochondria, it is recommended to isolate the mitochondria from fresh tissue or cells.

Add 5–50 μL of the samples into duplicate wells. Bring samples to a final volume of 50 μL using SCS Assay Buffer.

Note: For unknown samples, it is suggested to test several sample dilutions to ensure the readings are within the linear range of the standard curve.

For samples exhibiting significant background, especially background caused by NADH in the sample, include a Sample Blank for each sample by omitting the SCS Substrate Mix. The Sample Blank readings can then be subtracted from the sample readings.

For a positive control (optional), add 1–10 μL of the SCS Positive Control solution to the desired wells. Adjust the final volume to 50 μL with the SCS Assay Buffer.

Assay Reaction

1. Set up the Reaction Mixes according to the scheme in Table 1. 50 μL of Reaction Mix is required for each reaction (well).

Table 1.
Reaction Mixes

| Reagent | Standards, Controls, and Samples | Sample Blank |
|-------------------|----------------------------------|------------------|
| SCS Assay Buffer | 44 μL | 46 μL |
| SCS Substrate Mix | 2 μL | – |
| SCS Enzyme Mix | 2 μL | 2 μL |
| SCS Developer | 2 μL | 2 μL |

2. Add 50 μL of the appropriate Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting.
3. Measure the absorbance (A_{450}) in a microplate reader in kinetic mode for 10–30 minutes at $25\text{ }^{\circ}\text{C}$. Protect the plate from light during the incubation. It is recommended to take absorbance readings every minute.
Note: Incubation time depends on the activity of SCS in the samples.
4. Continue taking measurements until the value of the most active sample is greater than the value of the highest standard (12.5 nmole/well). At this time the most active sample is near or exceeds the end of the linear range of the standard curve.

Note: The NADH Standards can be read at the end of the incubation time.

Results

Calculations

Correct for the background by subtracting the measurement obtained for the 0 (blank) NADH Standard from that of the standards, controls, and samples. Background values can be significant and must be subtracted from all readings. Use the values obtained from the appropriate NADH Standards to plot a standard curve.

Note: A new standard curve must be set up each time the assay is run.

Plot the absorbance (A_{450}) for each well versus time.

Choose two time points (T1 and T2) in the linear range of the plot and determine the A_{450} at each time (ABS1 and ABS2).

Note: It is essential that ABS1 and ABS2 fall within the linear range of the standard curve.

Calculate the change in absorbance from T1 to T2 for the samples.

$$\Delta\text{ABS} = \text{ABS2} - \text{ABS1}$$

Subtract the Sample Blank ΔABS value from the Sample ΔABS reading to obtain the corrected measurement. Using the corrected measurement, determine the amount of NADH (nmole/well) generated by the SCS assay between T1 and T2 (S_a).

Succinyl-CoA Synthetase activity

$$\text{SCS Activity} = \frac{S_a}{(\text{Reaction Time}) \times S_v}$$

where:

S_a = Amount of NADH (nmole) generated in unknown sample well between T1 and T2 from standard curve

Reaction Time = T2 – T1 (minutes)

S_v = sample volume (μL) added to well

SCS activity is reported as nmole/min/ μL = milliunit/ μL .

One unit of succinyl-CoA synthetase is the amount of enzyme that generates 1.0 μmole of NADH per minute at pH 7.4 at 25 °C.

Sample Calculation

Amount of NADH (S_a) = 5.84 nmole
(from standard curve)

(T1) = 3 minutes

(T2) = 32 minutes

Sample volume (S_v) = 50 μL

SCS activity in sample well:

$$\text{nmole/min}/\mu\text{L} = \frac{5.84 \text{ nmole/well}}{(32 \text{ min} - 3 \text{ min}) \times 50 \mu\text{L/well}} = 0.00403$$

(milliunits/ μL)

References

1. Lamperti, C. et al., A novel homozygous mutation in SUCLA2 gene identified by exome sequencing. *Mol. Genet. Metab.*, **107**, 403–408 (2012).
2. Elpeleg, O. et al., Deficiency of the ADP-forming succinyl-CoA synthase activity is associated with encephalomyopathy and mitochondrial DNA depletion. *Am. J. Hum. Genet.*, **76**, 1081–1086 (2005).

Troubleshooting Guide

| Problem | Possible Cause | Suggested Solution |
|--|---|--|
| Assay not working | Cold assay buffer | Assay Buffer must be at room temperature |
| | Omission of step in procedure | Refer and follow Technical Bulletin precisely |
| | Plate reader at incorrect wavelength | Check filter settings of instrument |
| | Type of 96 well plate used | For colorimetric assays, use clear plates |
| Samples with erratic readings | Samples prepared in different buffer | Use the Assay Buffer provided or refer to Technical Bulletin for instructions |
| | Cell/Tissue culture samples were incompletely homogenized | Repeat the sample homogenization, increasing the length and extent of homogenization step. |
| | Samples used after multiple freeze-thaw cycles | Aliquot and freeze samples if needed to use multiple times |
| | Presence of interfering substance in the sample | If possible, dilute sample further |
| | Use of old or inappropriately stored samples | Use fresh samples and store correctly until use |
| Lower/higher readings in samples and standards | Improperly thawed components | Thaw all components completely and mix gently before use |
| | Use of expired kit or improperly stored reagents | Check the expiration date and store the components appropriately |
| | Allowing the reagents to sit for extended times on ice | Prepare fresh Reaction Mixes before each use |
| | Incorrect incubation times or temperatures | Refer to Technical Bulletin and verify correct incubation times and temperatures |
| | Incorrect volumes used | Use calibrated pipettes and aliquot correctly |
| Non-linear standard curve | Use of partially thawed components | Thaw and resuspend all components before preparing the Reaction Mixes |
| | Pipetting errors in preparation of standards | Avoid pipetting small volumes |
| | Pipetting errors in the Reaction Mix | Prepare Reaction Mixes whenever possible |
| | Air bubbles formed in well | Pipette gently against the wall of the plate well |
| | Standard stock is at incorrect concentration | Refer to the standard dilution instructions in the Technical Bulletin |
| | Calculation errors | Recheck calculations after referring to Technical Bulletin |
| | Substituting reagents from older kits/lots | Use fresh components from the same kit |
| Unanticipated results | Samples measured at incorrect wavelength | Check the equipment and filter settings |
| | Samples contain interfering substances | If possible, dilute sample further |
| | Sample readings above/below the linear range | Concentrate or dilute samples so readings are in the linear range |

GD,KVG,LS,MAM 10/15-1