

## ab287844 – Mitochondrial Complex III Activity Assay Kit

For in vitro quantitative determination of the activity of complex III in isolated mitochondria.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab287844>

### Storage and Stability

An unopened kit can be stored at -20°C for 6 months.

### Materials Supplied

Item	Quantity	Storage Condition
Complex III Assay Buffer	25 mL	-20°C
Oxidized Cytochrome c	4 X 1 vials	-20°C
Antimycin A	200 µL	-20°C
DTT IV	1 mL	-20°C
96-Well Half Area Plate	1 unit	-20°C

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Multi-well spectrophotometer capable of reading absorbance in kinetic mode
- DMSO (anhydrous)
- Deionized water

### Reagent Preparation

- Before using the kit, spin the tubes prior to opening.

Complex III Assay Buffer: Store at 4°C. Bring to RT before use.

Oxidized Cytochrome c: Reconstitute one vial at a time with 175 µL Complex III Assay Buffer to obtain 2 mM solution. Once reconstituted, one vial is enough for 25 reactions. Centrifuge briefly after mixing. Store at -20°C. The reconstituted vial stored at -20°C should be stable for at least one month.

Antimycin A: Aliquot and store at -20°C.

DTT IV: Thaw before use.

**Δ Note:** Antimycin A and DTT IV are stable for at least 3 months at -20°C.

### Sample Preparation

Isolate mitochondria from cultured cells or tissue using preferred procedure or kit.

- Estimate the protein concentration of isolated mitochondrial samples using Bradford assay.
- Isolated mitochondria should be stored at -80°C unless being used for the assay immediately. Avoid freeze-thaw cycles.
- Different dilutions of the mitochondrial sample should be tested to make sure that the activity falls in the linear range of the assay.
- Dilutions should be prepared in Complex III Assay Buffer immediately before performing the assay.

**Δ Note:** Mitochondria should be placed on ice during the assay.

Version 4b, Last updated Wednesday, June 18, 2025

### Assay Protocol

#### Reduced Cytochrome c Standard Curve

1. Prepare 750 mM DTT IV solution by mixing 1 M DTT IV solution and Assay Buffer in a ratio 3:1.
2. Add 0, 2, 4, 6, 8 and 10 µL of 2.0 mM Oxidized Cytochrome c standard into a series of wells in the provided 96-Well Half Area Plate to generate 0, 4, 8, 12, 16, and 20 nmol/well of Oxidized Cytochrome c.
3. Adjust the volume to 23 µL/well with Complex III Assay Buffer.
4. Add 2 µL of the prepared 750 mM DTT IV solution in each well in order to completely reduce Oxidized Cytochrome c.
5. Incubate at RT for 10 min on a plate shaker set to 200 rpm.
6. Measure the absorbance at 550 nm (end point).

#### Reaction Mix

- Mix enough reagents for the number of assays to be performed. For each well, prepare reaction mix containing:

Item	Background Control Well	Sample Mix Well	Sample Mix with Antimycin A
Assay Buffer	17 µL	15 µL	15 µL
Antimycin A	-	-	2 µL
DMSO	2 µL	2 µL	-

- Add the reaction mix to wells of the clear bottom 96-well half area plate which is provided with the kit.

#### Mitochondrial Sample

1. Add 1 to 2 µL mitochondrial samples (1.5 to 10 µg protein) to wells containing "Sample Mix" and "Sample Mix with Antimycin A". Mix well.
2. Add 6 µL Oxidized Cytochrome c (substrate) and read immediately.

**Δ Note:** Have the 96-Well Half Area Plate/plate reader ready at 550 nm on kinetic mode.

#### Measurement

- Immediately start recording absorbance at 550 nm, at 30 second intervals for up to 10 minutes at room temperature.

## Calculation

- Complex III specific activity may be calculated by comparing sample OD values (after subtraction of background control) to the reduced Oxidized Cytochrome c standard curve.
- Calculate the concentration of reduced Oxidized Cytochrome c, at time t1 and t2 by reading off the standard curve.
- Calculate  $\Delta$  [cytochrome c] between time t1 and t2. Apply the following equation to obtain activity of complex III:

$$\text{Complex III specific activity} = \frac{\Delta c}{\Delta t \times p} \times D \text{ (Units}/\mu\text{g)}$$

Where:

- $\Delta C$  is the change in reduced cytochrome c concentration during  $\Delta t$
  - $\Delta t$  is the time duration between t1 and t2, equal to: t2 – t1 (min)
  - p is the mitochondrial protein sample ( $\mu\text{g}$ )
  - D is the dilution factor
- 
- Use the following equation to calculate net activity:
- $$\text{Net Complex III Activity in sample} =$$
- $$\text{Activity in reaction without Antimycin A} - \text{Activity in reaction with Antimycin A}$$
- 
- Unit definition: One unit of Complex III is the amount of enzyme that will cause reduction of 1.0  $\mu\text{mol}$  of cytochrome c per min at pH 7.4 at room temperature.

## Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)