

**ab178779**

# **Yeast Mitochondria Isolation Kit**

## Instructions for Use

For the isolation of mitochondria from *P. pastoris* and other species of yeast

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

# Table of Contents

---

Table of Contents	1
1. Overview	2
2. Protocol Summary	3
3. Kits Components	4
4. Storage and Stability	5
5. Materials Required, Not Supplied	5
6. Reagents Preparation	6
7. Assay Protocol	7

# 1. Overview

---

Mitochondria are semi-autonomous organelles which functions in aging process, apoptosis, anti-HIV drugs, and cancers. Mitochondrial DNA (mtDNA) has a very high mutation rate and the mutations on mtDNA appear to be related to certain diseases such as diabetes, Alzheimer's disease, and muscle disorders. Isolation and quantification of mtDNA are often required to study the relationships between the diseases and mtDNA.

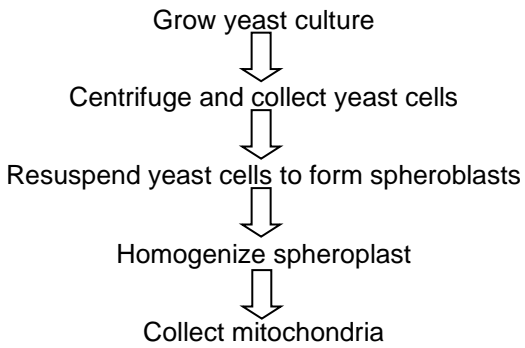
Abcam's Yeast Mitochondria Isolation Kit (ab178779) provides a convenient tool for isolating mitochondria from a variety of yeast species, and it has been specifically tested on *Pichia pastoris* and *Saccharomyces cerevisiae*.

Isolated mitochondria can be used for studying mitochondrial respiratory chain, assembly of complexes, apoptosis, mtDNA and mtRNA, and for proein profiling. Isolated mitochondria can be used directly in Western blot and ELISA.

ab178779 is only sufficient for 25 samples.

## 2. Protocol Summary

---



### 3. Kits Components

---

<b>Item</b>	<b>Quantity</b>
Buffer A	50 mL
Buffer B	50 mL
DTT IV	1 mL
Yeast Homogenization Buffer	50 mL
Lysis Enzyme Mix	220 $\mu$ L
Dilution Buffer I	10 mL
Protease Inhibitor Cocktail I	1 vial

PLEASE NOTE: DTT IV was previously labelled as 1 M DTT, and Yeast Homogenization Buffer as Homogenization Buffer, and Protease Inhibitor Cocktail I as Protease Inhibitor Cocktail (lyophilized), and Dilution Buffer I as Storage Buffer. The composition has not changed.

## **4. Storage and Stability**

---

Upon arrival, store the kit at -20°C and protect from light. Please read the entire protocol before performing the assay. Avoid repeated freeze/thaw cycles.

Briefly centrifuge all small vials prior to opening.

## **5. Materials Required, Not Supplied**

---

- Ultrapure water or MilliQ
- DMSO
- Appropriate media to grow yeast cells
- Glass douncer
- Spectrophotometer capable of reading absorbance
- Centrifuge with cooling option

## 6. Reagents Preparation

---

### 1. Dilution Buffer I:

Ready to use as supplied. Keep on ice while in ice. Store at -20°C or 4°C.

### 2. Protease Inhibitor Cocktail I (PIC):

Resuspend Protease Inhibitor Cocktail I in 250 µL of DMSO. Store at -20°C.

### 3. Lysis Enzyme Mix:

Aliquot Lysis Enzyme Mix and store at -20°C. Avoid repeated freeze/thaw.

### 4. DTT IV:

Ready to use as supplied. Store at -20°C.

### 5. Buffer A:

Store at -20°C or 4°C. Warm to room temperature before use. Prior to use, add DTT IV to a final concentration of 10mM to buffer A. Buffer A + DTT IV solution has to be used fresh and cannot be stored for future uses.

### 6. Buffer B:

Store at -20°C or 4°C. Warm to room temperature before use.

Prior to use, add Lysis Enzyme 5 $\mu$ L/mL of buffer B. Buffer + Lysis Enzyme solution has to be used fresh and cannot be stored for future uses.

#### **7. Yeast Homogenization Buffer:**

Store at -20°C or 4°C. Prior to use, add Protease Inhibitor Cocktail I to 1:1000 final dilution. Yeast Homogenization Buffer + Protease Inhibitor Cocktail I solution has to be used fresh and cannot be stored for future uses.

## **7. Assay Protocol**

---

This procedure is for small-scale isolation (10 – 20 mL) for OD~20. For a large scale preparation (OD~200), calculate the volumes of the reagents accordingly.

#### **1. Yeast Culture:**

- a)** Grow yeast cells in appropriate media overnight at 30°C, shaking at 200 rpm. For temperature sensitive mutants, use the desired temperature.
- b)** When cells are into log phase, determine the OD of the culture at 600 nm. Multiply the OD with the total volume of the culture (mL) to calculate the total OD.

*NOTE: To isolate mitochondria in respiring state, grow yeast cells under aerobic condition using non-fermentable media (e.g. Ethanol*

*or Glycerol as carbon source). However, yeast cells will grow very slowly under these conditions with a thicker cell wall.*

## **2. Mitochondrial Isolation:**

- a)** Centrifuge the yeast culture at 3,000 rpm for 5 min and discard the supernatant. Wash the cells by resuspending in 2 volumes of ultrapure water.
- b)** Resuspend the cell pellet in 1 mL of Buffer A containing 10 mM fresh DTT IV and incubate for 10 min at 30°C with gentle shaking. Centrifuge at 1,500 g for 5 min and discard the supernatant.
- c)** Resuspend the cell pellet in 1 mL of Buffer B. Aliquot 10 µL suspension in a separate glass tube (Control).
- d)** Add 5 µL Lysis Enzyme Mix to the remaining cell suspension and incubate for 10 – 15 min at 30°C in a shaking incubator. Aliquot 10 µL of suspension again in another glass tube.

*NOTE: To check the formation of efficient spheroplast (cell from which the cell wall has been almost completely removed), add 990 µL of water to 10 µL aliquots of step 2c and 2d (Control & with Lysis Enzyme Mix). Measure OD at 600 nm. Incubation should continue until the OD of the sample is decreased 30 – 40% after adding Lysis Enzyme Mix compared to Control.*

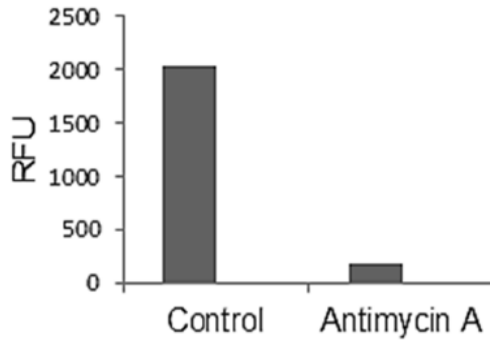
- e)** After efficient spheroplast formation, centrifuge at 1,500 g for 5 min and discard the supernatant. From this step onwards, keep tubes on ice.

- f) Resuspend the cell pellet in 1 mL of Yeast Homogenization Buffer + Protease Inhibitor Cocktail I. Transfer the suspension to a glass douncer and stroke 10 – 15 times on ice. Centrifuge at 600 g for 5 min at 4°C and collect the supernatant in a separate tube. The supernatant contains the mitochondria.
- g) Centrifuge the supernatant containing mitochondria again at 600 g for 5 min at 4°C and collect the supernatant.
- h) Centrifuge the supernatant at 12,000 g for 10 min at 4°C. Carefully discard the supernatant without touching the pellet.
- i) Resuspend the pellet in ~ 50 µL of Dilution Buffer I. Determine the protein concentration and adjust the desired protein concentration by Dilution Buffer I accordingly.

### **3. Isolated Mitochondria Storage:**

The storage conditions will vary based on the application the samples will be used for:

- Intact mitochondria: resuspend mitochondria in Dilution Buffer I and snap freeze in liquid nitrogen. Transfer frozen mitochondria to storage at -80°C.
- Mitochondria for Gel loading: mitochondria can be stored in Lysis Buffer with detergent or SDS-PAGE loading dye.



**Figure 1:** Mitochondrial integrity test – Purified mitochondria were analyzed for intactness by using JC-1 dye, which tests the electrochemical proton gradient ( $\Delta\Psi$ ) of the inner mitochondrial membrane. The intact purified mitochondria show aggregation of JC-1 dye whose signal can be measured at Ex/Em = 530/590 nm. Treatment with antimycin A (100  $\mu$ M), an inhibitor of the electron transport chain, dissipates the mitochondrial membrane potential resulting in reduced fluorescence signal.

## **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)