

# Mitochondria Isolation Kit for Cultured Cells

89874

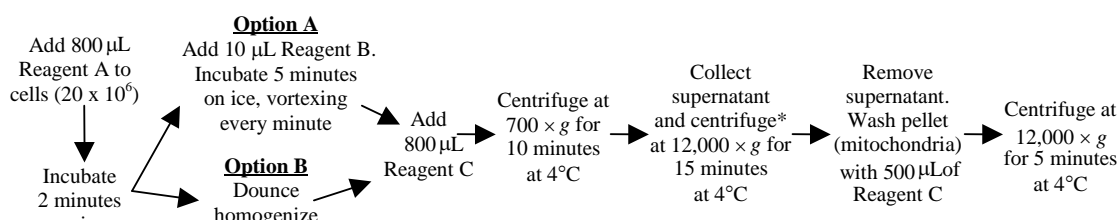
1477.7

Number	Description
89874	<b>Mitochondria Isolation Kit for Cultured Cells</b> , contains sufficient reagents to isolate intact mitochondria from 50 pellets of cultured mammalian cells containing $2 \times 10^7$ cells each
	<b>Kit Contents:</b>
	<b>Mitochondria Isolation Reagent A</b> , 50mL
	<b>Mitochondria Isolation Reagent B</b> , 500 $\mu$ L
	<b>Mitochondria Isolation Reagent C</b> , 65mL

**Storage:** Upon receipt store kit at 4°C. Product is shipped at ambient temperature.

## Introduction

The Thermo Scientific™ Mitochondria Isolation Kit for Cultured Cells enables isolation of intact mitochondria from cultured mammalian cells. The kit offers two options for the separation of mitochondria from cytosolic components (Figure 1). One option is a unique protocol utilizing a reagent-based method allowing multiple samples to be processed concurrently. The second option uses traditional Dounce homogenization and provides approximately two-fold more mitochondria, based on protein analysis. Both protocols rely on differential centrifugation to separate the mitochondrial and cytosolic fractions with a bench-top microcentrifuge and are completed in approximately 40 minutes (post-cell harvest). Each procedure has been optimized for maximum yield of mitochondria with minimal damage to integrity. Once isolated, the mitochondria may be used for a number of downstream applications, including apoptosis, signal transduction and metabolic studies.



\*To obtain a more purified preparation of mitochondria, this step may be performed at 3,000  $\times$  g.

**Figure 1.** Procedure summary of the Pierce Mitochondria Isolation Kit.

## Important Product Information

- This kit allows two options for mitochondria isolation: a reagent-based method or Dounce homogenization. The kit contains sufficient reagents for a **total of 50 applications** using either method exclusively or a combination of both.
- Up to six samples may be processed concurrently using the reagent-based isolation method, whereas only one sample may be processed at a time with the Dounce isolation method. Because of the time delay in processing multiple samples with the Dounce protocol, samples processed initially may yield a higher quantity of mitochondria than cell pellets incubated on ice for subsequent isolation.

- The protocols presented prepare a crude mitochondrial fraction that includes lysosomal and peroxisomal contaminants. To obtain a more purified fraction of heavy mitochondria, centrifuge the post-nuclear supernatant at  $3000 \times g$  instead of at  $12,000 \times g$ . Western blot analysis of purified versus crude mitochondrial fractions prepared with this kit result in a  $> 50\%$  reduction in cathepsin S (lysosomal protein) and PMP70 (peroxisomal protein) in the purified fraction.

## Additional Materials Required

- Variable-speed bench-top microcentrifuge (refrigerated)
- 2.0mL microcentrifuge tubes
- Vortex mixer
- Protease inhibitors, EDTA-free such as Thermo Scientific™ Halt™ Protease Inhibitor Cocktail, EDTA-Free (100X), 1mL (Product No. 87785)
- Dounce tissue grinder, such as 2mL Kontes or Wheaton Dounce Tissue Grinder, if using Option B

## Option A: Isolation of Mitochondria using Reagent-based Method

### Notes:

- Immediately before use, add protease inhibitors to Reagent A and Reagent C; only add inhibitors to the reagent amount being used for the procedure and not to the stock solutions.
- Process up to six samples concurrently.
- Required speed of vortex changes during the protocol.

1. Pellet  $2 \times 10^7$  cells by centrifuging harvested cell suspension in a 2.0mL microcentrifuge tube at  $\sim 850 \times g$  for 2 minutes. Carefully remove and discard the supernatant.
2. Add 800 $\mu$ L of Mitochondria Isolation Reagent A. Vortex at **medium** speed for 5 seconds and incubate tube on ice for exactly 2 minutes.

**Note:** Do not exceed the 2 minute incubation.

3. Add 10 $\mu$ L of Mitochondria Isolation Reagent B. Vortex at **maximum** speed for 5 seconds.
4. Incubate tube on ice for 5 minutes, vortexing at **maximum** speed every minute.
5. Add 800 $\mu$ L of Mitochondria Isolation Reagent C. Invert tube several times to mix (**do not vortex**).
6. Centrifuge tube at  $700 \times g$  for 10 minutes at  $4^\circ\text{C}$ .
7. Transfer the supernatant to a new, 2.0mL tube and centrifuge at  $12,000 \times g$  for 15 minutes at  $4^\circ\text{C}$ .

**Note:** To obtain a more purified fraction of mitochondria, with  $> 50\%$  reduction of lysosomal and peroxisomal contaminants, centrifuge at  $3000 \times g$  for 15 minutes.

8. Transfer the supernatant (cytosol fraction) to a new tube. The pellet contains the isolated mitochondria.
9. Add 500 $\mu$ L Mitochondria Isolation Reagent C to the pellet, and centrifuge at  $12,000 \times g$  for 5 minutes. Discard the supernatant.
10. Maintain the mitochondrial pellet on ice before downstream processing. Freezing and thawing may compromise mitochondria integrity.

## Option B: Isolation of Mitochondria using Dounce Homogenization

### Notes:

- Immediately before use, add protease inhibitors to Reagent A and Reagent C; only add inhibitors to the reagent amount being used for the procedure and not to the stock solutions.
- Process one sample at a time.
- Pre-chill Dounce Tissue Grinder on ice before use.

1. Pellet  $2 \times 10^7$  cells by centrifuging harvested cell suspension in a 2.0mL microcentrifuge tube at  $\sim 850 \times g$  for 2 minutes. Carefully remove and discard the supernatant.
2. Add 800 $\mu$ L of Mitochondria Isolation Reagent A. Vortex at **medium** speed for 5 seconds and incubate tube on ice for exactly 2 minutes.  
**Note:** Do not exceed the 2 minute incubation.
3. Transfer cell suspension to Dounce Tissue Grinder.
4. Homogenize cells on ice. Perform enough strokes to effectively lyse the cells.  
**Note:** See Additional Information Section for the number of strokes required for  $> 80\%$  lysis of a select group of cell types.  
**Note:** To check the cell lysis efficiency, spot 5 $\mu$ L of cell lysate onto a glass slide, add coverslip and view under a microscope. Compare with 5 $\mu$ L of the non-lysed cells.
5. Return lysed cells to original tube and add 800 $\mu$ L of Mitochondria Isolation Reagent C.
6. Rinse Dounce Tissue Grinder with 200 $\mu$ L of Mitochondria Isolation Reagent A and add to tube containing the sample in step B.5.
7. Invert tube several times to mix (**do not vortex**).
8. Centrifuge tube at  $700 \times g$  for 10 minutes at  $4^\circ\text{C}$ .
9. Transfer the supernatant to a new, 2.0mL tube and centrifuge at  $12,000 \times g$  for 15 minutes at  $4^\circ\text{C}$ .  
**Note:** To obtain a more purified fraction of mitochondria, with  $> 50\%$  reduction of lysosomal and peroxisomal contaminants, centrifuge at  $3000 \times g$  for 15 minutes.
10. Transfer the supernatant (cytosol fraction) to a new tube. The pellet contains the isolated mitochondria.
11. Add 500 $\mu$ L Mitochondria Isolation Reagent C to the pellet, and centrifuge at  $12,000 \times g$  for 5 minutes. Discard the supernatant.
12. Maintain the mitochondrial pellet on ice before downstream processing. Freezing and thawing may compromise mitochondria integrity.

## Additional Information

### A. Cell Lysis

The number of Dounce homogenization strokes necessary for optimal cell lysis will vary depending upon cell line. The number of strokes required for  $> 80\%$  lysis of a select group of cell types is indicated in Table 1, which may be used as a guide for other cell lines.

**Table 1.** Number of strokes required to achieve  $\sim 80\%$  lysis efficiency using Dounce homogenization.\*

<u>Cell Line</u>	<u>Number of strokes</u>
C6	40
HeLa	60
NIH 3T3	80

\* Lysis efficiency was determined by visual estimation using a microscope.

### B. Mitochondria Lysis

For analysis by Western blotting or gel electrophoresis, boil mitochondrial pellet with SDS-PAGE sample buffer and apply to the gel. For protein analysis using the Thermo Scientific™ BCA Protein Assay Kit (Product No. 23225), mitochondria may be lysed with 2% CHAPS in Tris-buffered saline (TBS; e.g., 25mM Tris, 0.15M NaCl; pH 7.2; Product No. 28379) as described below:

1. Add 100 $\mu$ L of 2% CHAPS in TBS to the mitochondrial pellet and vortex for 1 minute.
2. Centrifuge mitochondria at high speed for 2 minutes. The supernatant contains soluble mitochondrial protein that can be analyzed by BCA Protein Assay.

---

## Related Thermo Scientific Products

<b>89801</b>	<b>Mitochondria Isolation Kit for Tissue</b>
<b>23225</b>	<b>BCA Protein Assay Kit</b>
<b>34080</b>	<b>SuperSignal™ West Pico Chemiluminescent Substrate, 500mL</b>
<b>87785</b>	<b>Halt™ Protease Inhibitor Cocktail, EDTA-Free (100X), 1mL</b>
<b>78833</b>	<b>NE-PER™ Nuclear and Cytoplasmic Extraction Kit</b>
<b>89842</b>	<b>Mem-PER™ Plus Membrane Protein Extraction Kit</b>

## References for Apoptosis

Green, D.R. and Reed, J.C. (1998). Mitochondria and apoptosis. *Science* **281**:1309-12.

Newmeyer, D.D. and Ferguson-Miller, S. (2003). Mitochondria: Releasing power for life and unleashing the machineries of death. *Cell* **112**:481-90.

## References for Mitochondrial Proteome

Lescuyer, P., *et al.* (2003). Progress in the definition of a reference human mitochondrial proteome. *Proteomics* **3**:157-67.

Taylor, S.W., *et al.* (2003). Characterization of the human heart mitochondrial proteome. *Nat Biotechnol* **21**:281-5.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.