

Protocol for Total RNA Isolation from Cells

Description

The following protocol is for total RNA isolation from cells using the Total RNA Purification Kit (Catalog #79040). For complete instructions, refer to the Technical Manual (Document #10000005434).

Directions

1. Prepare cell lysate from adherent cells or cells in suspension, as indicated below.

Adherent Cells

- a) Aspirate cell culture medium. Add ice-cold sterile D-PBS as indicated in Table 1. Aspirate D-PBS.

Table 1. Recommended Volumes of PBS, RNA Lysis Buffer + TG, and 100% Isopropanol for Various Cultureware

Cultureware	Volume of D-PBS	Volume of RNA Lysis Buffer + TG	Volume of 100% Isopropanol
T-25 cm ² flask	5 mL	500 µL	170 µL
6-well plate	2 mL	250 µL	85 µL
24-well plate	500 µL	100 µL	35 µL
48-well plate	250 µL	100 µL	35 µL
96-well plate	100 µL	100 µL	35 µL

- b) Add RNA Lysis Buffer + TG as indicated in Table 1. Gently rock the plate or flask to completely cover the adherent cells with buffer. Pipette the lysate up and down over the cultureware surface 7 - 10 times.
- c) Collect the lysate and transfer to a new microcentrifuge tube.
- d) Add 100% isopropanol as indicated in Table 1. Mix by vortexing for 5 seconds.
- e) Proceed to step 2 for RNA isolation.

Cells in Suspension

- a) In a sterile centrifuge tube, centrifuge cell suspension at 300 x g for 5 minutes. Remove and discard supernatant.

- b) Add ice-cold, sterile D-PBS to wash cells. Centrifuge at 300 x g for 5 minutes. Remove as much supernatant as possible and discard.
- c) Add RNA Lysis Buffer + TG as indicated in Table 2. Mix well by pipetting up and down 7 - 10 times, or by vortexing. For > 2 x 10⁶ cells, pass the lysate through a 20-gauge needle 4 - 5 times to shear the genomic DNA.
- d) Add 100% isopropanol as indicated in Table 2. Mix by vortexing for 5 seconds.
- e) Proceed to step 2 for RNA isolation.

Table 2. Recommended Volumes of RNA Lysis Buffer + TG and Isopropanol per Cell Input Range

Cell Input Range	Volume of RNA Lysis Buffer + TG	Volume of 100% Isopropanol
1 x 10 ² to 5 x 10 ⁵	100 µL	35 µL
> 5 x 10 ⁵ to 2 x 10 ⁶	250 µL	85 µL
> 2 x 10 ⁶ to 5 x 10 ⁶	500 µL	170 µL

2. Insert minicolumn into Collection Tube.
3. Transfer the lysate to the minicolumn assembly.
4. Centrifuge at 12,000 - 14,000 x g for 30 seconds. Discard the liquid in the Collection Tube and reinsert minicolumn into Collection Tube.
5. Add 500 µL RNA Wash Buffer to the minicolumn. Centrifuge at 12,000 - 14,000 x g for 30 seconds. Discard the liquid in the Collection Tube and reinsert minicolumn into Collection Tube.
6. Prepare DNase I Incubation Mix by combining the reagents as indicated in Table 3, per sample, **in the order listed**.

Table 3. Preparation of DNase I Incubation Mix

Component	Volume Per Sample
DNA Digestion Buffer	24 µL
MnCl ₂	3 µL
DNase I Solution	3 µL

7. Mix by gently pipetting up and down; **do not vortex**. Store on ice.
8. Add 30 µL of fresh DNase I Incubation Mix directly to the minicolumn membrane. Incubate at room temperature (15 - 25°C) for 15 minutes.
9. Add 200 µL of Column Wash Buffer (with ethanol added) to the minicolumn. Centrifuge at 12,000 - 14,000 x g for 15 seconds.
10. Add 500 µL of RNA Wash Buffer (with ethanol added) to the minicolumn. Centrifuge at 12,000 - 14,000 x g for 30 seconds. Remove the minicolumn and transfer to a new Collection Tube. Discard the Collection Tube containing wash buffer.
11. Add 300 µL of RNA Wash Buffer to the minicolumn. Centrifuge at high speed for 2 minutes.
12. Transfer minicolumn to an Elution Tube. Add nuclease-free water to the minicolumn membrane as indicated in Table 4. Incubate at room temperature for 1 - 2 minutes. Centrifuge at 12,000 - 14,000 x g for 1 minute.
13. Remove and discard minicolumn. Store purified RNA at -80°C.

Table 4. Recommended Volume of Nuclease-Free Water per Cell Input Range

Cell Input Range	Volume of Nuclease-Free Water
1 x 10 ² to 5 x 10 ⁵	15 µL
> 5 x 10 ⁵ to 2 x 10 ⁶	30 µL
> 2 x 10 ⁶ to 5 x 10 ⁶	50 µL

Copyright © 2019 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, and Scientists Helping Scientists are trademarks of STEMCELL Technologies Canada Inc. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

STEMCELL TECHNOLOGIES INC.'S QUALITY MANAGEMENT SYSTEM IS CERTIFIED TO ISO 13485. PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES UNLESS OTHERWISE STATED.

Protocol for Total RNA Isolation

