

RNA Isolation and Purification Procedure from Cell Samples

Use the following protocol to lyse cultured cells from suspension. Use 1×10^2 to 1×10^6 cells per purification.

1. Collect cells by centrifugation at $300 \times g$ for 5 minutes. Wash the cell pellet with ice-cold, sterile 1X PBS. Centrifuge at $300 \times g$ for 5 minutes to collect the cells. Discard the supernatant.
2. Add 200 μ l of LBA + TG Buffer to the washed cell pellet. Mix well by vortexing and/or pipetting.
Note: Following lysis, pipet 7–10 times to shear the DNA using a P200 or P1000 pipettor.
3. Add 130 μ l of RDB to each homogenate and vortex for 10 seconds. Centrifuge for 2 minutes at $12,000 \times g$. Carefully transfer the cleared homogenate to a clean 1.5ml tube.
4. Add 400 μ l of 100% isopropanol to each cleared homogenate. Mix by vortexing.
5. Transfer the homogenate to a ReliaPrep™ Minicolumn. Centrifuge at $12,000 \times g$ for 30 seconds.
6. Remove the column and discard the liquid. Place the column back into the Collection Tube.
7. Transfer the remaining homogenate to the same column used in Step 5. Centrifuge at $12,000 \times g$ for 30 seconds.
8. Remove the column and discard the liquid. Place the column back into the Collection Tube.
9. Add 500 μ l of RWA to each column. Centrifuge at $12,000 \times g$ for 30 seconds.
10. Remove the column and discard the liquid. Place the column back into the Collection Tube.
11. Add 500 μ l of RWA to each column. Centrifuge at $12,000 \times g$ for 2 minutes. Carefully transfer the column to a 1.5ml Elution Tube.
12. Add 40 μ l of Nuclease-Free Water to each column. Centrifuge at $12,000 \times g$ for 1 minute.
13. Transfer 5 μ l of DNase 10X Buffer and 5 μ l of DNase I to eluate.
14. Incubate for 5 minutes at room temperature (20–25°C).
15. Add 150 μ l of LBA + TG Buffer to the DNase treatment tube.
16. Add 300 μ l of 95% ethanol to the mixture and vortex for 10 seconds. Transfer 500 μ l of this mixture to a new column. Centrifuge at $12,000 \times g$ for 30 seconds.
17. Remove the column and discard the liquid. Place the column back into the Collection Tube and repeat Steps 9–11.
18. Add 15 μ l of Nuclease-Free Water to each column (see Table 1). Centrifuge at $12,000 \times g$ for 1 minute.

Table 1. Recommended RNA Elution Volumes per Number of Cells.

Cell Input Range	Nuclease-Free Water
1×10^2 to 5×10^5	15 μ l
$>5 \times 10^5$ to 1×10^6	30 μ l

Note: RNA concentration may increase with lower elution volumes; however, the total yield of RNA may decrease when elution volumes are between 10–15 μ l. If maximum recovery of RNA is essential, we recommend a second elution into a second sterile tube with an additional 15 μ l of Nuclease-Free Water followed by centrifugation at $12,000 \times g$ for 1 minute.

ReliaPrep™ miRNA Cell and Tissue Miniprep System

Instructions for Use of Products Z6210, Z6211, and Z6212.



Quick Protocol

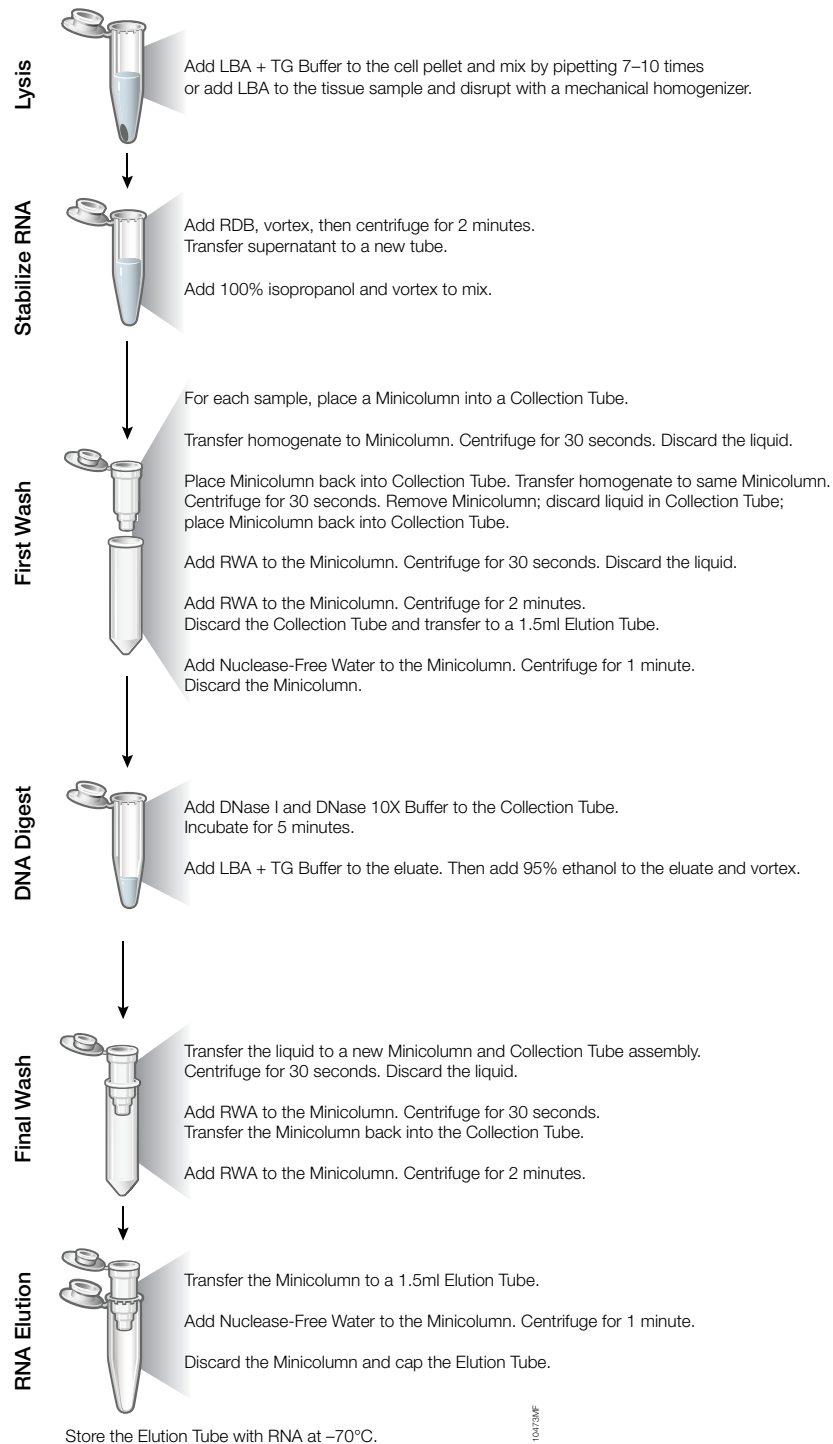


Figure 1. Schematic diagram of the ReliaPrep™ miRNA Cell and Tissue Miniprep System.

Additional protocol information is in Technical Manual #TM469, available online at: www.promega.com