

# ReliaPrep™ RNA Miniprep System for Cells

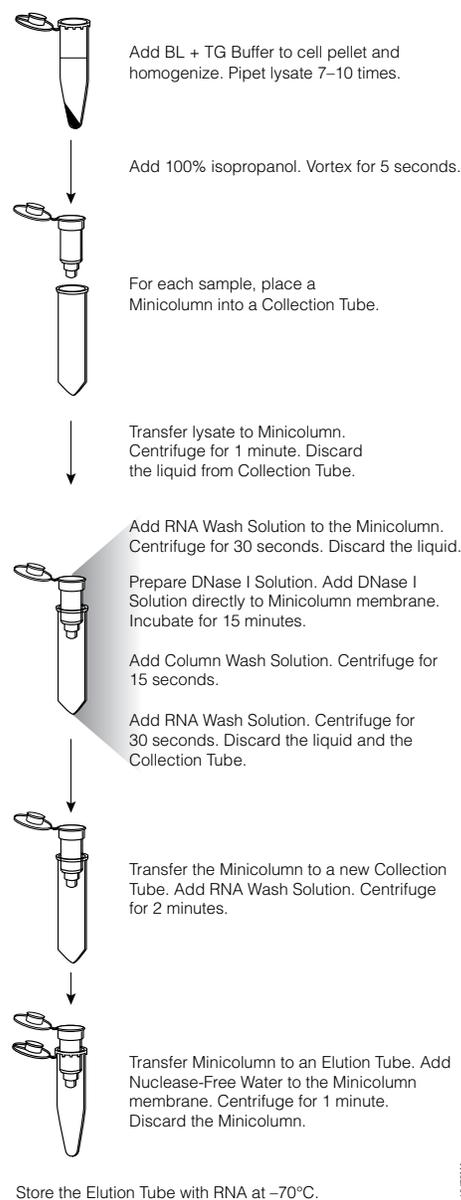
Instructions for Use of Products **Z6014** and **Z6015**.

## Protocol

- Before beginning the ReliaPrep™ RNA Miniprep for Cells protocol, see **Section 4.A., Preparing Solutions** in the *ReliaPrep™ RNA Miniprep System Technical Manual #TM774*. Prepare the four required solutions immediately prior to use, for best results.
- To collect and lyse adherent cells, use the protocol in Section 4.B, TM774, and then proceed to Step 7. For suspension cells, proceed to Step 3.
- Collect cells in a sterile centrifuge tube 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. Carefully discard the supernatant.
- Add BL + TG Buffer to the washed cell pellet (see the table below). If frozen cell pellets are used as starting material, add BL + TG Buffer to the frozen pellets before thawing.

Number of Cells	BL + TG Buffer	100% Isopropanol
$1 \times 10^2$ to $5 \times 10^5$	100µl	35µl
$>5 \times 10^5$ to $2 \times 10^6$	250µl	85µl
$>2 \times 10^6$ to $5 \times 10^6$	500µl	170µl

- Disperse the cell pellet and mix well by vortexing and/or pipetting.  
**Note:** After adding BL + TG Buffer, pipet 7–10 times to shear the DNA. For  $>2 \times 10^6$  cells, pass the lysate through a 20-gauge needle four to five times to shear the genomic DNA.
- Add isopropanol as recommended in the table above. Mix by vortexing 5 seconds.
- Wearing gloves, unpack one Minicolumn, two Collection Tubes and one Elution Tube for each sample. Label each tube and Minicolumn. Place one Minicolumn into a Collection Tube for each sample.
- Transfer lysate from Step 7 to a Minicolumn in a Collection Tube. Centrifuge at  $12,000\text{--}14,000 \times g$  for 30 seconds at  $20^\circ\text{--}25^\circ\text{C}$ .
- Remove the ReliaPrep™ Minicolumn and discard liquid in the Collection Tube. Replace the Minicolumn in the Collection Tube. Add **500µl of RNA Wash Solution** to the Minicolumn. Centrifuge at  $12,000\text{--}14,000 \times g$  for 30 seconds. Empty the Collection Tube.



# ReliaPrep™ RNA Miniprep System for Cells

Instructions for Use of Products **Z6014** and **Z6015**.

11. Prepare **DNase I incubation mix** by combining the following amounts of reagent, per sample, in the order listed:

Solution	Volume	× Number of Preps	= Total
Yellow Core Buffer	24µl		
MnCl <sub>2</sub> , 0.09M	3µl		
DNase I	3µl		

Mix by gently pipetting; **do not vortex**. The volumes listed above make enough DNase I mix for a single sample. Multiply this amount by the number of samples to calculate the amount of DNase I mix to prepare.

12. Apply **30µl of DNase I incubation mix** to the Minicolumn membrane. Incubate for 15 minutes at 20–25°C.
13. Add **200µl of Column Wash Solution** (with ethanol added) to the Minicolumn. Centrifuge at 12,000–14,000 × *g* for 15 seconds.
14. Add **500µl of RNA Wash Solution** (with ethanol added). Centrifuge at 12,000–14,000 × *g* for 30 seconds. Discard the wash solutions and the Collection Tube.
15. Place the ReliaPrep™ Minicolumn into a new Collection Tube. Add **300µl of RNA Wash Solution** and centrifuge at high speed for 2 minutes.
16. Transfer the ReliaPrep™ Minicolumn from the Collection Tube to an Elution Tube. Add **Nuclease-Free Water** to the Minicolumn membrane as recommended in the table below. Place the Minicolumn and Elution Tube into a centrifuge with the Elution Tube lid facing to the outside. Centrifuge at 12,000–14,000 × *g* for 1 minute.

Number of Cells	Nuclease-Free Water
1 × 10 <sup>2</sup> to 5 × 10 <sup>5</sup>	15µl
>5 × 10 <sup>5</sup> to 2 × 10 <sup>6</sup>	30µl
>2 × 10 <sup>6</sup> to 5 × 10 <sup>6</sup>	50µl

17. Discard the Minicolumn. Cap the Elution Tube containing the purified RNA and store at –70°C.

ReliaPrep is a trademark of Promega Corporation.

Additional protocol information is in Technical Manual #TM774, available online at: [www.promega.com](http://www.promega.com)