

TECHNICAL MANUAL

ReliaPrep™ RNA Clean-Up and Concentration System

Instructions for Use of Products
Z1071, Z1072 and Z1073



ReliaPrep™ RNA Clean-Up and Concentration System

All technical literature is available at: www.promega.com/protocols/
 Visit the web site to verify that you are using the most current version of this Technical Manual.
 E-mail Promega Technical Services if you have questions on use of this system: techserv@promega.com

1.	Description.....	1
2.	Product Components and Storage Conditions	2
3.	General Considerations	3
	3.A. Processing Capacity	3
	3.B. Downstream Applications	3
	3.C. Creating a Ribonuclease-Free Environment	3
4.	RNA Purification and/or Concentration	4
	4.A. Preparing Solutions	4
	4.B. Purifying and/or Concentrating RNA Samples	5
5.	Determining RNA Yield and Purity	6
6.	Troubleshooting.....	7
7.	Related Products.....	8

1. Description

The purity and concentration of RNA prepared from conventional isolation protocols (organic extraction or column-based procedures) is often affected by carryover of purification reagents like guanidine salts and ethanol. Furthermore, the requirements of some purification kits for large elution volumes can leave the user with RNA too dilute for some downstream applications. Additionally, a simpler and faster protocol for enzymatic RNA reactions (reverse transcription or post-DNase treatment) is also desired. The ReliaPrep™ RNA Clean-Up and Concentration System was designed to address these needs while maintaining good recovery of RNA samples and ensuring their usefulness in downstream applications such as RT-qPCR, Northern blot analysis or RNA sequencing. An added benefit is that dilute samples or sample pools up to 300µl can be processed over a single column and eluted in 15µl of water or TE buffer. You can process up to 12 samples in less than 10 minutes.



2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT #
ReliaPrep™ RNA Clean-Up and Concentration System	250 preps	Z1073

For Research Use. Each system contains sufficient reagents for 250 RNA purifications. Includes:

- 5 packs Reliaprep™ Minicolumns (50/pack)
- 1 pack Collection Tubes (250/pack)
- 5 packs Elution Tubes (50/pack)
- 2 × 20ml Membrane Binding Solution
- 24ml Column Wash Solution (CWE)
- 58.8ml RNA Wash Solution (RWA)
- 25ml Nuclease-Free Water

PRODUCT	SIZE	CAT #
ReliaPrep™ RNA Clean-Up and Concentration System	50 preps	Z1072

For Research Use. Each system contains sufficient reagents for 50 RNA purifications. Includes:

- 1 pack Reliaprep™ Minicolumns (50/pack)
- 1 pack Collection Tubes (50/pack)
- 2 packs Elution Tubes (25/pack)
- 20ml Membrane Binding Solution
- 5ml Column Wash Solution (CWE)
- 11.8ml RNA Wash Solution (RWA)
- 1.25ml Nuclease-Free Water

PRODUCT	SIZE	CAT #
ReliaPrep™ RNA Clean-Up and Concentration System	10 preps	Z1071

For Research Use. Each system contains sufficient reagents for 10 RNA purifications. Includes:

- 1 pack Reliaprep™ Minicolumns (10/pack)
- 1 pack Collection Tubes (10/pack)
- 2 packs Elution Tubes (5/pack)
- 4ml Membrane Binding Solution
- 1ml Column Wash Solution (CWE)
- 11.8ml RNA Wash Solution (RWA)
- 1.25ml Nuclease-Free Water

Storage Conditions: All components can be stored at 15–30°C.

3. General Considerations

3.A. Processing Capacity

Each column is capable of binding as much as 80µg of RNA. However, if concentrating or cleaning up RNA preparations of greater than 10µg, we recommend eluting twice or increasing elution volume.

3.B. Downstream Applications

RNA purified using the ReliaPrep™ RNA Clean-Up and Concentration System is suitable for use in many molecular biology applications such as RT-qPCR, Northern blotting or next-generation sequencing.

For all downstream applications, continue to protect your samples from RNases by wearing gloves and using solutions and tubes that are RNase-free. Using an RNase inhibitor such as Recombinant RNasin® Ribonuclease Inhibitor (Cat.# N2511) can ensure protection from nucleases that might be introduced during downstream processing.

3.C. Creating a Ribonuclease-Free Environment

Ribonucleases are extremely difficult to inactivate. Avoid inadvertently introducing RNases into your RNA during or after the purification. This is especially important if the RNA sample was difficult to obtain or is irreplaceable. The following notes may help you to prevent accidental RNase contamination of your sample:

1. Two of the most common sources of RNase contamination are the user's hands or bacteria and molds present on airborne dust particles. To prevent contamination from these sources, use sterile technique when handling the reagents supplied with the kit. Wear gloves at all times and use eye protection.
2. Whenever possible, sterile disposable plasticware should be used for handling RNA. These materials are generally RNase-free and thus do not require pretreatment to inactivate RNases. Sterile elution tubes are provided.
3. Reagents contain irritants and should be properly disposed of as hazardous waste. Do not use products containing bleach for sanitation.
4. Change pipette tips frequently to prevent sample cross-contamination.



4. RNA Purification and/or Concentration

For best results, do not stop the procedure once samples are loaded onto the column. After elution, store samples at -70°C . If desired, RNA can be eluted in TE buffer (not provided).

Materials to be Supplied by the User

- 100% isopropanol, RNase-free
- 95–100% ethanol, RNase-free
- 1.5ml microcentrifuge tubes
- microcentrifuge capable of maintaining $10,000 \times g$ (14,000rpm)

4.A. Preparing Solutions

Solution	Preparation	Notes
Column Wash Solution (CWE)	250 prep size (Part# A254C): Add 36ml of 95–100% ethanol to the bottle containing 24ml of concentrate. 50 prep size (Cat.# A254B): Add 7.5ml of 95–100% ethanol to the bottle containing 5ml of concentrate. 10 prep size (Part# A254A): Add 1.5ml of 95–100% ethanol to the bottle containing 1ml of concentrate.	After adding ethanol, mark on the bottle label that this step has been performed. The reagent is stable at $15-30^{\circ}\text{C}$ when tightly capped.
RNA Wash Solution (RWA)	250 prep size (Part# Z3091): Add 100ml of 95–100% ethanol to the bottle containing 58.8ml of concentrate. 50 prep size (Part# Z309B): Add 20ml of 95–100% ethanol to the bottle containing 11.8ml of concentrate. 10 prep size (Part# Z309B): Add 20ml of 95–100% ethanol to the bottle containing 11.8ml of concentrate.	After adding ethanol, mark on the bottle label that this step has been performed. The reagent is stable at $15-30^{\circ}\text{C}$ when tightly capped.
Membrane Binding Solution	Check that no precipitate is present in the bottle.	If there is precipitate, warm the contents in a 37°C waterbath, swirling until the precipitate redissolves.

4.B. Purifying and/or Concentrating RNA Samples

Notes:

1. This protocol requires 100% isopropanol (not provided).
2. Perform all centrifugation steps at $10,000 \times g$ (14,000rpm).

Dilute RNA (μ l)	Membrane Binding Solution (μ l)	100% Isopropanol (μ l)	Total Load (μ l)
25	12.5	37.5	75
50	25	75	150
100	50	150	300
150	75	225	450
200	100	300	600
250	125	375	750
300	150	450	900

Note: You can process up to 300 μ l of dilute RNA at one time in a single minicolumn. If a dilute RNA sample exceeds 300 μ l, process the sample in multiple tubes in the manner described above and load the multiple centrifugations onto the same column. The wash protocol remains unchanged.

1. Pipet 25–300 μ l of dilute RNA into a 1.5ml microcentrifuge tube.
2. Add 1/2 volume of Membrane Binding Solution and vortex 5 seconds.
3. Add 1 1/2 volumes of 100% isopropanol.
4. Load RNA sample onto a ReliaPrep™ Minicolumn seated in a Collection Tube and centrifuge for 30 seconds.
Note: This mixture is viscous so slowly transfer the sample to the column. Save the sample tubes after transfer to check that no significant volume remains.
5. Remove column, and discard the contents of the Collection Tube. Reseat the column into the same Collection Tube.
6. Add 200 μ l of Column Wash Solution (CWE) and centrifuge for 15 seconds. Remove column, and discard the contents of the Collection Tube. Reseat the column into the same Collection Tube.
7. Wash with 300 μ l of RNA Wash Solution (RWA) and centrifuge for 15 seconds. Repeat wash with 300 μ l of RNA Wash Solution (RWA) and centrifuge again.
8. Remove column and discard the contents of the collection tube. Reseat the column into the same collection tube, centrifuge for 1 minute to dry the column and then transfer column to an Elution Tube.
9. Pipet 15 μ l of Nuclease-Free Water or TE buffer (not provided) to the center of the column matrix and centrifuge for 30 seconds.
Note: Touch the pipette tip to the column bed surface before dispensing Nuclease-Free Water or TE buffer to completely wet the column matrix. The color should change from light to dark tan.
10. For maximum recovery, repeat elution with an additional 15 μ l of Nuclease-Free Water or TE buffer.

5. Determining RNA Yield and Purity

The ReliaPrep™ RNA Clean-Up and Concentration System can be used to purify and concentrate RNA sourced from previous purification schemes or simply concentrate RNA preparations from dilute solutions. The total RNA yield may be determined spectrophotometrically at 260nm, where 1 absorbance unit (A_{260}) equals approximately 40µg of single-stranded RNA/ml. The purity also may be estimated by spectrophotometry from the relative absorbances at 230nm, 260nm and 280nm (i.e., A_{260}/A_{280} and A_{260}/A_{230} ratios). If overall expected yield is less than 1µg, spectrophotometric analysis will not yield accurate results due to the lack of absorbance sensitivity for low-concentration nucleic acids. Alternatively, low RNA yield can be determined with good accuracy using the QuantiFluor® RNA System (Cat. # E3310).

RNA purified using the ReliaPrep™ RNA Clean-Up and Concentration System is substantially free of contaminating protein. Pure RNA should exhibit an A_{260}/A_{280} ratio of 2.0. However, variations between individual starting materials and skill in performing the procedure may reduce this purity ratio to the 1.7–2.1 range. This is not necessarily problematic for downstream applications. RNA will usually exhibit an A_{260}/A_{230} ratio of 1.8–2.2. A low A_{260}/A_{230} ratio can indicate guanidine contamination, which may interfere with performance or accuracy of results in downstream applications. A low A_{260}/A_{230} ratio may also indicate ethanol contamination. Be sure to follow the centrifugation recommendations in the protocol.

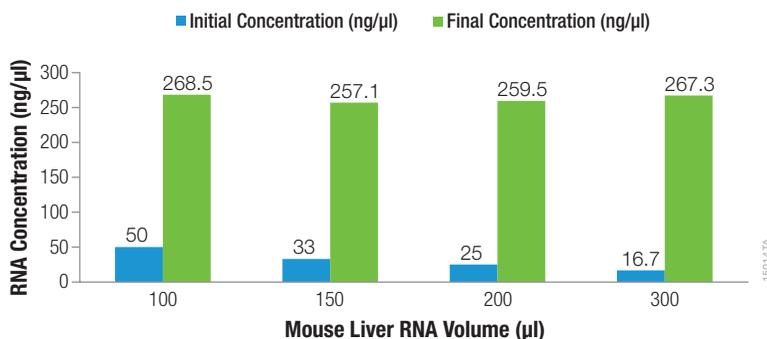


Figure 1. Recovery and concentration of RNA from dilute solution using the ReliaPrep™ RNA Clean-Up and Concentration System. Five micrograms of mouse liver RNA was suspended in 100, 150, 200 or 300µl of water, repurified using the standard protocol and eluted in 15µl of Nuclease-Free Water. Results shown in Figure 1 demonstrate equivalent yield regardless of target RNA concentration. Average RNA yield was 79.8%.

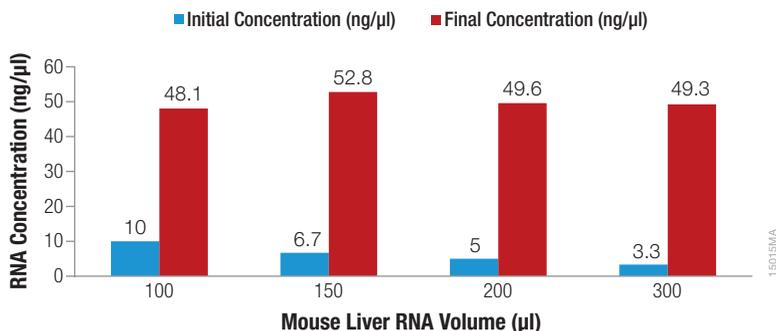


Figure 2. Recovery and concentration of RNA from dilute solution using the ReliaPrep™ RNA Clean-Up and Concentration System. One microgram of mouse liver RNA was suspended in 100, 150, 200 or 300μl of water, repurified using the standard protocol and eluted in 15μl of Nuclease-Free Water. Results shown in Figure 2 demonstrate equivalent yield regardless of target RNA concentration. Average RNA yield was 74.9%.

6. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com. E-mail: techserv@promega.com

Symptoms

Low A_{260}/A_{280} ratios

Possible Causes and Comments

Low RNA concentration. Spectrophotometric methods may not be accurate when expected RNA concentration drops below approximately 15ng/μl. As the signal approaches the limit of detection for the instrument, background noise may lead to inaccurate readings.

Low A_{260}/A_{230} ratios

Spectrophotometric methods may not be accurate when expected RNA concentration drops below approximately 15ng/μl. As the signal approaches the limit of detection for the instrument, background noise may lead to inaccurate readings.

Low A_{260}/A_{230} ratios are typically due to guanidine thiocyanate contamination. Be certain to complete all wash steps in the protocol.

Low A_{260} value

Low RNA yield. Spectrophotometric methods may not be accurate when expected RNA concentration drops below approximately 15ng/μl. As the signal approaches the limit of detection for the instrument, background noise may lead to inaccurate readings.



6. Troubleshooting (continued)

Symptoms	Possible Causes and Comments
Low A ₂₆₀ value (continued)	<p>The correct reagent volume may not have been added during the protocol. The volumes used for the kit reagents are critical to the success of the procedure.</p> <p>Ethanol may not have been added to the Column Wash Solution or RNA Wash Solution. Prepare solutions as instructed in Section 4.A.</p>
Calculated yield lower than expected	RNA may be retained on the minicolumn if expected yield is >10µg. Elute a second time with 15µl of Nuclease-Free Water.
RNA degradation	<p>RNase was introduced during handling. Carefully handle kit reagent bottles to avoid contamination. Use the provided disposable plasticware when manipulating and storing RNA. Wear gloves at all times.</p> <p>RNA was degraded during sample preparation. Work quickly when preparing samples, and store eluates at -70°C.</p>

7. Related Products

Product	Size	Cat.#
PureYield™ Plasmid Miniprep System	100 preps	A1223
	250 preps	A1222
PureYield™ Plasmid Midiprep System	25 preps	A2492
	100 preps	A2495
	300 preps	A2496
PureYield™ Plasmid Maxiprep System	10 preps	A2392
	25 preps	A2393
pGEM®-T Vector Systems	20 reactions	A3600
pGEM®-T Vector Systems + JM109 Competent Cells	20 reactions	A3610
pGEM®-T Easy Vector Systems	20 reactions	A1360
pGEM®-T Easy Vector Systems + JM109 Competent Cells	20 reactions	A1380
LigaFast™ Rapid DNA Ligation System	30 reactions	M8221
	150 reactions	M8225
QuantiFluor® RNA System	1ml	E3310

Note: You can use the CloneWeaver™ Workflow Builder to design a cloning workflow for your construct:
www.promega.com/resources/tools/cloneweaver/

© 2018 Promega Corporation. All Rights Reserved.

pGEM, QuantiFluor and RNasin are registered trademarks of Promega Corporation. CloneWeaver, LigaFast, PureYield and ReliaPrep are trademarks of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.