# Maxwell<sup>®</sup> 16 LEV simplyRNA Cells Kit and Maxwell<sup>®</sup> 16 LEV simplyRNA Tissue Kit

Instructions for Use of Products AS1270 and AS1280

Promega

Note: To run the simplyRNA protocol, you must have Maxwell<sup>®</sup> 16 firmware version ≥4.95 (Cat.# AS2000) or ≥1.50 (Cat.# AS3000) installed on your instrument, and you must use the Maxwell<sup>®</sup> 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070). Using the original LEV magnetic rod will result in low yields.





# Maxwell<sup>®</sup> 16 LEV simplyRNA Cells Kit and Maxwell<sup>®</sup> 16 LEV simplyRNA Tissue Kit

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.	Before You Begin	3 4 4
4.	Purification of Total RNA from Tissue Samples	6
5.	Purification of Total RNA from Cell Culture	7
6.	Instrument Run: AS2000 and AS3000 Instruments   6.A. Setup for AS2000 Maxwell® 16 Instruments   6.B. Setup for AS3000 Maxwell® 16 MDx Instruments	8 8 10
7.	Troubleshooting	. 12
8.	Appendix   8.A. Creating a Ribonuclease-Free Environment   8.B. Downstream Applications   8.C. References   8.D. Related Products	14 14 14 14 15
	8.E. Summary of Changes	. 16
	• •	

## 1. Description

The Maxwell<sup>®</sup> 16 LEV simplyRNA Cells Kit<sup>(a)</sup> (Cat.# AS1270) and Maxwell<sup>®</sup> 16 LEV simplyRNA Tissue Kit<sup>(a)</sup> (Cat.# AS1280) are used with the Maxwell<sup>®</sup> 16 Instrument (Cat.# AS2000 or AS3000) configured with the Maxwell<sup>®</sup> 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070) to purify RNA from fresh or frozen mammalian cells or fresh or frozen mammalian tissue samples, respectively. This RNA purification procedure is a simple method with minimal lysate handling before automated purification on the Maxwell<sup>®</sup> 16 Instrument. The low elution volume is used to generate concentrated high-quality RNA suitable for use in downstream applications such as quantitative RT-PCR. The kits provide the reagents for processing the samples and use prefilled cartridges for purification, maximizing simplicity and convenience. The Maxwell<sup>®</sup> 16 Instrument can process from 1 to 16 samples in under an hour.



## 2. Product Components and Storage Conditions

PRO	DUCT		SIZE	CAT.#
Maxwell <sup>®</sup> 16 LEV simplyRNA Cells Kit		LEV simplyRNA Cells Kit	48 preps	AS1270
For	Laborate	ory Use. Sufficient for 48 automated isolations from cell samples. Inclu	ıdes:	
•	48	Maxwell® 16 LEV Cartridges (MCE)		
•	30ml	Homogenization Solution		
•	20ml	Lysis Buffer		
•	1 vial	DNase I (lyophilized)		
•	900µl	1-Thioglycerol		
•	50µl	Blue Dye		
•	25ml	Nuclease-Free Water		
•	50	LEV Plungers		
•	50	Elution Tubes, 0.5ml		

PRODUCT	SIZE	CAT.#
Maxwell <sup>*</sup> 16 LEV simplyRNA Tissue Kit	48 preps	AS1280

For Laboratory Use. Sufficient for 48 automated isolations from tissue samples. Includes:

- 48 Maxwell<sup>®</sup> 16 LEV Cartridges (MCE)
- 30ml Homogenization Solution
- 20ml Lysis Buffer
- 2 vials DNase I (lyophilized)
- 900µl 1-Thioglycerol
- 50µl Blue Dye
- 25ml Nuclease-Free Water
- 50 LEV Plungers
- 50 Elution Tubes, 0.5ml

**Storage Conditions:** Upon receipt, remove 1-Thioglycerol and store at  $2-10^{\circ}$ C. Store the remaining kit components at room temperature (15–30°C). 1-Thioglycerol also can be stored at room temperature (15–30°C), where it is stable for up to 9 months.

**Safety Information:** The reagent cartridges contain ethanol, which is flammable. 1-Thioglycerol is toxic. Guanidine thiocyanate and guanidine hydrochloride (which are components of the Homogenization Solution and Lysis Buffer) are harmful and irritants. The Lysis Buffer also has a possible risk of harm to an unborn child. Wear gloves and follow standard safety procedures while working with these substances.



The Maxwell<sup>®</sup> 16 reagent cartridges are designed to be used with potentially infectious substances. Wear appropriate protection (e.g., gloves and goggles) when handling infectious substances. Adhere to your institutional guidelines for the handling and disposal of all infectious substances used with this system.

**Note:** Bleach reacts with guanidine thiocyanate and should not be added to any sample waste containing the Homogenization Solution.

# 3. Before You Begin

# 3.A. Maxwell® 16 Instrument Hardware and Firmware Setup

To use the Maxwell<sup>®</sup> 16 LEV simplyRNA Tissue Kit or Cells Kit, the Maxwell<sup>®</sup> 16 Instrument must be configured with LEV hardware. If your Maxwell<sup>®</sup> 16 Instrument contains standard elution volume (SEV) hardware or the original LEV Magnetic Rod, **it must be reconfigured using the Maxwell<sup>®</sup> 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070**; Figure 1). Using the original LEV magnetic rod will result in low yields. Reconfiguring the instrument is simple. Refer to the *Maxwell<sup>®</sup> 16 Instrument Operating Manual* specific for your instrument for directions.



# Figure 1. Maxwell 16<sup>®</sup> High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070).

Failure to change the Maxwell<sup>®</sup> 16 Instrument hardware to the LEV configuration could result in instrument damage. Using the standard elution volume (SEV) hardware configuration with LEV-configured reagent products will cause damage to the instrument.

**Important:** To run the simplyRNA protocol, you must have Maxwell<sup>®</sup> 16 firmware version  $\geq$ 4.95 (Cat.# AS2000) or  $\geq$ 1.50 (Cat.# AS3000) installed on your instrument.

**Capacity:** The Maxwell<sup>®</sup> 16 simplyRNA Tissue Kit can process up to 20mg of most tissues. Higher amounts of some tissues may result in higher yields (e.g., heart). The Maxwell<sup>®</sup> 16 simplyRNA Cells Kit can process up to 5 million cells.

## 3.B. Preparation of Solutions

## **Homogenization Solution**

To prepare a working solution, add 20 $\mu$ l of 1-Thioglycerol per milliliter of Homogenization Solution. 1-Thioglycerol is viscous, so careful pipetting is required for accurate measurement. Alternatively, add 600 $\mu$ l of 1-Thioglycerol to the 30ml bottle of Homogenization Solution. A volume of 200 $\mu$ l of 1-Thioglycerol/Homogenization Solution is needed for each sample. Before use, chill the 1-Thioglycerol/Homogenization Solution on ice or at 2–10°C.

Note: Store the 1-Thioglycerol/Homogenization Solution at 2–10°C, where it is stable for up to 30 days.

## DNase I

Add 275 $\mu$ l of Nuclease-Free Water to the vial of lyophilized DNase I. Invert to rinse DNase off the underside of the cap and swirl gently to mix; do not vortex. Add 5 $\mu$ l of Blue Dye to the reconstituted DNase I as a visual aid for pipetting. Dispense the DNase I solution into single-use aliquots in nuclease-free tubes. Store reconstituted DNase I at  $-20^{\circ}$ C. Do not freeze-thaw reconstituted DNase I more than three times.

## **3.C. Cartridge Preparation**

To maintain an RNase-free environment during processing, change gloves before handling cartridges, LEV Plungers and Elution Tubes. Place the cartridges to be used in the Maxwell<sup>®</sup> 16 LEV Cartridge Rack (Cat.# AS1251). Place each cartridge in the rack with the label side facing away from the Elution Tubes. Press down on the cartridge to snap it into position. Carefully peel back the seal so that all plastic comes off the top of the cartridge. Ensure that all sealing tape and any residual adhesive are removed before placing cartridges in the instrument.

Note: If you are processing fewer than 16 samples, center the cartridges on the platform.

- 1. Place an LEV Plunger in well #8 of each cartridge. Well #8 is the well closest to the Elution Tube.
- 2. Place 0.5ml Elution Tubes in the front of the Maxwell<sup>®</sup> 16 LEV Cartridge Rack. Add 50µl of Nuclease-Free Water to the bottom of each Elution Tube. For a more concentrated eluate, as little as 30µl of nuclease-free water may be added to the elution tube, but the total amount of RNA recovered may be reduced.

#### Notes:

- 1. If Nuclease-Free Water is on the side of the tube, the elution may be suboptimal.
- 2. Use only the 0.5ml Elution Tubes provided in the kit; other tubes may not work with the Maxwell® 16 Instrument.

For setup of tissue samples, see Section 4. For setup of cell culture samples, see Section 5.



Figure 2. The Elution Tubes are placed in the front of the Maxwell® 16 LEV Cartridge Rack, and 50µl of Nuclease-Free Water is dispensed into each tube.



Figure 3. The LEV plunger is placed in well #8 of the cartridge (the well closest to the Elution Tube), and lysates are placed into well #1 of the cartridge.



## 4. Purification of Total RNA from Tissue Samples

#### Materials to Be Supplied By the User

- small tissue homogenizer (e.g., Tissue-Tearor™ homogenizer, PRO Scientific or any homogenizer capable of handling small volumes)
- vortex mixer
- tube for homogenization
- RNase-free, sterile, aerosol-resistant pipette tips
- **optional:** heat block or water bath set to 70°C
- 1. Working as quickly as possible, homogenize the tissue sample in the chilled 1-Thioglycerol/Homogenization Solution (Section 3.B) until no visible tissue fragments remain. Homogenize an additional 15–30 seconds for complete homogenization. If foaming occurs, let sample settle on ice. Extra Homogenization Solution is provided, but only 200µl of homogenate can be processed per cartridge. The final volume of the homogenate to be added to the cartridge should be 200µl. Add additional homogenization solution as needed to bring samples to a final volume of 200µl.

**Note:** Samples may be stored frozen at  $-80^{\circ}$ C after homogenization for later processing. Thaw homogenates on ice or at  $2-10^{\circ}$ C to avoid RNA degradation.

2. **Optional:** RNA yield from larger amounts of liver may be increased by heating homogenates at 70°C for 2 minutes, then allowing homogenates to cool (approximately 1 minute) before proceeding to Step 3. This is recommended for 10mg or more of liver tissue.

**Note:** If the heat step is used, the purified RNA may be partially denatured, and may migrate differently on native gels. Denaturing gels are recommended if the heating step is used.

- 3. Shortly before processing samples on the Maxwell<sup>®</sup> 16 Instrument, add 200µl of Lysis Buffer (Part #MC501C) to 200µl of homogenate. Vortex vigorously for 15 seconds to mix. Transfer all 400µl of lysate to well 1 of the Maxwell<sup>®</sup> 16 LEV Cartridge (MCE). Well #1 is closest to the cartridge label and farthest from the elution tube.
- 4. Add 5µl of DNase (Section 3.B) to well 4 (yellow reagent). When using more than 5mg of tissues with high DNA content (e.g., liver or spleen), add 10µl of DNase to well 4. After the blue DNase I solution is added, the reagent in well 4 will be green.

Proceed to Section 6 for instructions on loading samples on the instrument and beginning the automated purification.



## 5. Purification of Total RNA from Cell Culture

#### Materials to Be Supplied By the User

- centrifuge
- vortex mixer
- RNase-free, sterile, aerosol-resistant pipette tips
- 1. Trypsinize adherent cells following normal protocols.
- 2. Pellet cells at low speed (e.g.,  $300 \times g$  for 3 minutes).
- 3. Remove medium.
- 4. Add 200µl of chilled 1-Thioglycerol/Homogenization Solution (Section 3.B) to the cell pellet and vortex until pellet is dispersed and cells appear lysed. A pipette may be used to disperse pellets before vortexing. Alternatively cells can be homogenized. Store lysed cells on ice if there is a delay before processing.
- 5. Shortly before processing samples on the Maxwell<sup>®</sup> 16 Instrument, add 200µl of Lysis Buffer (Part# MC501C) to 200µl of lysed cells. Vortex vigorously for 15 seconds to mix. Transfer all 400µl of lysate to well 1 of the Maxwell<sup>®</sup> 16 LEV Cartridge (MCE). Well #1 is the closest to the cartridge label and farthest from the elution tube.
- 6. Add 5μl of DNase I solution to well 4 (yellow reagent). After adding the blue DNase, the solution in well 4 will be green.



#### 6. Instrument Run: AS2000 and AS3000 Instruments

#### 6.A. Setup for AS2000 Maxwell® 16 Instruments

Refer to the Maxwell® 16 Instrument Operating Manual #TM295 for more detailed information.

**Important:** To run the simplyRNA protocol, you must have Maxwell<sup>®</sup> 16 firmware version **4.95** or higher installed on your instrument, and you must use the Maxwell<sup>®</sup> 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070). Using the original LEV magnetic rod will result in low yields.

1. Turn on the Maxwell<sup>®</sup> 16 Instrument. The instrument will power up, display the firmware version number, proceed through a self-check and home all moving parts.

Note: Ensure that you are running firmware version 4.95 or higher.

- 2. Verify that the instrument settings indicate an "LEV" hardware configuration and "Rsch" operational mode setting.
- 3 Select "Run" on the Menu screen, and press the Run/Stop button to select the method.
- 4. Select "RNA", select "simplyRNA", then select "simplyRNA" once more on the Menu screen. Next select "OK" at the Verification screen.
- 5. Open the door when prompted. Press the Run/Stop button to extend the platform.



Warning: Pinch point hazard.

6. Transfer the Maxwell<sup>®</sup> 16 LEV Cartridge Rack containing prepared cartridges on the Maxwell<sup>®</sup> 16 Instrument platform. Ensure that the rack is placed in the Maxwell<sup>®</sup> 16 Instrument with the Elution Tubes closest to the door. The rack will only fit in the instrument in this orientation. If you have difficulty fitting the rack on the instrument platform, check that the rack is in the correct orientation. Ensure that the cartridge rack is level on the instrument platform.

Note: Hold the Maxwell® 16 LEV Cartridge Rack by the sides to avoid dislodging cartridges from the rack.

7. Verify that samples were added to well #1 of the cartridges, cartridges in the rack are loaded on the instrument, Elution Tubes are present with 50µl of Nuclease-Free Water and LEV Plungers are in well #8. Well #4 should be green to indicate that DNase was added.



Note: Failure to add DNase will result in DNA in the eluate.

8. Press the Run/Stop button. The platform will retract. Close the door.



Warning: Pinch point hazard.

9. The Maxwell<sup>®</sup> 16 Instrument will immediately begin the purification run. The screen will display the steps performed and the approximate time remaining in the run.

## Notes:

- 1. Pressing the Run/Stop button or opening the door will pause the run.
- 2. If the run is abandoned before completion, the instrument will wash the particles off the plungers and eject the plungers into well #8 of the cartridge. To continue processing the samples, rinse any particles off the plunger into the last well used. Discard the used plungers. Put new plungers into well #8, and start the run from the beginning.
- 10. When the automated purification run is complete, the LCD screen will display a message that the method has ended.

## End of Run

- 11. Follow the on-screen instructions at the end of the method to open door. Verify that the plungers are located in well #8 of the cartridge at the end of the run. If the plungers are not removed from the magnetic plunger bar, push them down by hand to remove them.
- 12. Press the Run/Stop button to extend the platform out of the instrument.
- 13. Remove the Maxwell<sup>®</sup> 16 LEV Cartridge Rack from the instrument. Remove Elution Tubes containing total RNA, and close the tubes.
- 14. If paramagnetic particles are present in the elution tubes, centrifuge at  $10,000 \times g$  for 2 minutes.

Alternatively, if desired, an additional particle capture step may be performed using the 0.5ml MagneSphere<sup>®</sup> Technology Magnetic Separation Stand (Cat.# Z5341) or Maxwell<sup>®</sup> 16 LEV Magnet (Cat.# AS1261). Transfer the supernatant to a clean tube (not provided). Avoid transferring paramagnetic particles.

15. Remove the cartridges and plungers from the Maxwell<sup>®</sup> 16 LEV Cartridge Rack, and discard following the recommended guidelines. Do not reuse reagent cartridges, LEV plungers or Elution Tubes.

## **Storing Eluted RNA**

If sample eluates are not processed immediately, the eluted RNA should be stored at  $-70^{\circ}$ C, or at  $-20^{\circ}$ C for up to 24 hours. Consult the protocol for your downstream application for specific storage and handling recommendations.

## 6.B. Setup for AS3000 Maxwell® 16 MDx Instruments

Refer to the Maxwell® 16 MDx Instrument Technical Manual #TM320 for detailed information.

**Important:** To run the simplyRNA protocol, you must have Maxwell<sup>®</sup> 16 firmware version ≥**1.50** (AS3000) installed on your instrument, and you must use the Maxwell<sup>®</sup> 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070). Using the original LEV magnetic rod will result in low yields.

- 1. Turn on the Maxwell<sup>®</sup> 16 MDx Instrument. The instrument will power up, display the firmware version number, proceed through a self-check and home all moving parts.
- 2. Verify that the Home screen indicates "LEV" and the LEV hardware is present. Press "Run" to continue.
- 3 Enter user and PIN if this option is enabled.
- 4. At the Protocols screen, select "RNA", then select "simplyRNA".
- 5. On the next screen, verify that the correct method and user were chosen. Select "Run/Stop" to continue.
- 6. Open the door when prompted on the screen, then select "Run/Stop".



Warning: Pinch point hazard.

- 7. Follow on-screen instructions for bar code reader input if this option is enabled.
- 8. Transfer the Maxwell<sup>®</sup> 16 LEV Cartridge Rack containing prepared cartridges on the Maxwell<sup>®</sup> 16 Instrument platform. Ensure that the rack is placed in the Maxwell<sup>®</sup> 16 Instrument with the Elution Tubes closest to the door. The rack will only fit in the instrument in this orientation. If you have difficulty fitting the rack on the instrument platform, check that the rack is in the correct orientation. Ensure that the cartridge rack is level on the instrument platform.

Note: Hold the Maxwell<sup>®</sup> 16 LEV Cartridge Rack by the sides to avoid dislodging cartridges from the rack.

9. Verify that samples were added to well #1 of the cartridges, cartridges in the rack are loaded on the instrument, Elution Tubes are present with 50µl of Nuclease-Free Water and LEV Plungers are in well #8. Well #4 should be green to indicate that DNase was added.



Note: Failure to add DNase will result in DNA in the eluate.

10. Press the Run/Stop button. The platform will retract. Close the door.



Warning: Pinch point hazard.

The Maxwell<sup>®</sup> 16 MDx Instrument will immediately begin the purification run. The screen will display the approximate time remaining in the run.

#### Notes:

- 1. Pressing the Run/Stop button or opening the door will pause the run.
- 2. If the run is abandoned before completion, the instrument will wash the particles off the plungers and eject the plungers into well #8 of the cartridge. To continue processing the samples, rinse any particles off the plunger into the last well used. Discard the used plungers. Put new plungers into well #8, and start the run from the beginning.
- 11. When the automated purification run is complete, follow instructions on the screen for data transfer. For detailed instructions, refer to the *Maxwell*<sup>®</sup> 16 MDx Instrument Technical Manual #TM320 and Maxwell<sup>®</sup> Sample Track Software Technical Manual #TM314.

## End of Run

- 12. Follow on-screen instructions at the end of the method to open door. Verify that plungers are located in well #8 of the cartridge at the end of the run. If plungers are not removed from the magnetic plunger bar, push them down gently by hand to remove them.
- 13. Press the Run/Stop button to extend the platform out of the instrument.
- 14. Remove the Maxwell® 16 LEV Cartridge Rack from the instrument. Remove Elution Tubes containing RNA, and close the tubes.
- 15. If paramagnetic particles are present in the elution tubes, centrifuge at  $10,000 \times g$  for 2 minutes.

Alternatively, if desired, an additional particle capture step may be performed using the 0.5ml MagneSphere<sup>®</sup> Technology Magnetic Separation Stand (Cat.# Z5341) or Maxwell<sup>®</sup> 16 LEV Magnet (Cat.# AS1261). Transfer the supernatant to a clean tube (not provided). Avoid transferring paramagnetic particles.

16. Remove the cartridges and plungers from the Maxwell<sup>®</sup> 16 LEV Cartridge Rack, and discard following the recommended guidelines. Do not reuse reagent cartridges, LEV plungers or Elution Tubes.

For the Maxwell® 16 MDx Instrument, ensure samples are removed before the UV light treatment to avoid damage to the nucleic acid.

## **Storing Eluted RNA**

If sample eluates are not processed immediately, the eluted RNA should be stored at  $-70^{\circ}$ C, or at  $-20^{\circ}$ C for up to 24 hours. Consult the protocol for your downstream application for specific storage and handling recommendations.

# 7. 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	Causes and Comments
Sample foams during homogenization	Some homogenizers will generate foam when samples are homogenized. Allow the foam to dissipate prior to pipetting. Homogenize for shorter periods of time until visible particles and tissue fragments are eliminated. Keep rotor submerged whenever the homogenizer is on.
	Sample was homogenized in the Lysis Buffer instead of the Homogenization Solution.
Homogenate is too viscous to pipet	The homogenate was too concentrated and became viscous while sitting on ice. Reduce the homogenate viscosity by increasing the amount of Homogenization Solution 1.5- to 2-fold, and briefly rehomogenize the sample. The maximum volume of homogenate that can be processed in a single Maxwell <sup>®</sup> Cartridge (MCE) is 200µl.
Low RNA yield, RNA degradation or	1-Thioglycerol was not added to the Homogenization Solution.
poor reproducibility between samples	Lysis Buffer was not added.
	Lysates were not mixed by vortexing long enough.
	Homogenization was incomplete. Incomplete homogenization of samples results in loss of RNA within the particulates and clumps of debris.
	RNA yield for liver may be improved by incubation at 70°C for 2 minutes.
	Samples were not properly prepared or stored. Samples must be flash frozen, stored in RNA <i>later</i> <sup>®</sup> reagent or immediately homogenized in Homogenization Solution to halt RNA degradation. Delays during sample collection may result in RNA degradation and lower yields. Freeze samples immediately, and store at -70°C if they cannot be processed immediately. <u>Homogenates should be stored at -70°C and thawed on ice</u> . Frozen lysate was heated to thaw. Thaw frozen lysates on ice or
	at 2–10°C.
	The wrong magnet bar was used. Use the High Strength LEV Magnetic Rod and Plunger Bar Adaptor (Cat.# SP1070).

12 Promega Corporation · 2800 Woods Hollow Road · Madison, WI 53711-5399 USA · Toll Free in USA 800-356-9526 · 608-274-4330 · Fax 608-277-2516 TM351 · Revised 8/17 www.promega.com

Symptoms	Causes and Comments		
Low RNA yield, RNA degradation or poor reproducibility between samples (continued)	Sample contains a low amount of RNA. The amount of RNA present in a sample depends on the metabolic state, stage of growth, type of sample and growth conditions. Sample types vary in the amount of total RNA.		
	RNase introduced by handling. Use sterile, disposable plastic- ware or baked glassware when handling RNA. Wear clean gloves at all times. RNases introduced during or after purification will degrade the RNA. See Section 8.A, Creating a Ribonuclease- Free Environment.		
	The wrong method was run with the Maxwell® 16 Instrument.		
DNA contamination seen when performing RT-PCR or PCR	DNase was not added to the correct well in the cartridge, or DNase not added at all. Check the color of the liquid in well 4. If the blue DNase Solution was added, the reagent in well 4 will be green, not yellow.		
	Too much sample was processed. Reduce the starting sample amount twofold.		
	Sample has an excessive amount of genomic DNA. Reduce the starting material or increase the amount of DNase added.		
	Possible cross-contamination. RT-PCR and PCR are extremely sensitive techniques. Use aerosol-resistant pipette tips. Set up reactions and analyze samples in separate locations.		
	Too much sample was used in RT-PCR. Reduce the total RNA input to 50–100ng in RT-PCR. Generally a rare message can be detected in 50ng of total RNA by RT-PCR.		
	The wrong method was run with the Maxwell® 16 Instrument.		
Purified total RNA appears cloudy	Total RNA purified from liver may contain glycogen. When stored at 4°C or frozen, the glycogen may form a precipitate, and the sample may appear cloudy. Warm the sample to 23–25°C, and vortex to dissolve the glycogen. Glycogen does not interfere in reactions that use nucleic acids as a substrate.		
In a gel, eluate floats out of the well when loading	Alcohol carryover in the eluate may cause it to float. Allow eluate to air dry, or speed vac before loading in a gel.		



## 8. Appendix

#### 8.A. Creating a Ribonuclease-Free Environment

Ribonucleases are extremely difficult to inactivate. Take care to avoid introducing RNase activity into your RNA samples during and after isolation. This is especially important if the starting material was difficult to obtain or is irreplaceable. The following notes may help prevent accidental RNase contamination of your samples.

- 1. Two of the most common sources of RNase contamination are the user's hands and bacteria or molds that may be present on airborne dust particles. To prevent contamination from these sources, use sterile technique when handling the reagents supplied with this system. Wear gloves at all times. Change gloves whenever ribonucleases may have been contacted.
- 2. Whenever possible, sterile, disposable plasticware should be used for handling RNA. These materials generally are RNase-free and do not require pretreatment to inactivate RNase.
- 3. Treat nonsterile glassware, plasticware and electrophoresis chambers before use to ensure that they are RNasefree. Bake glassware at 200°C overnight, and thoroughly rinse plasticware with 0.1N NaOH, 1mM EDTA, followed by RNase-free water. Commercially available RNase removal products also may be used, following the manufacturer's instructions.

**Note:** Electrophoresis chambers may be contaminated with ribonucleases, particularly RNase A, from analysis of DNA samples. Whenever possible, set aside a new or decontaminated apparatus for RNA analysis only.

4. Treat solutions not supplied with the system by adding diethyl pyrocarbonate (DEPC) to 0.1% in a fume hood. Incubate overnight with stirring at room temperature in the hood. Autoclave for 30 minutes to remove any trace of DEPC.

**Caution:** DEPC is a suspected carcinogen and should only be used in a chemical fume hood. DEPC reacts rapidly with amines and cannot be used to treat Tris buffers.

#### 8.B. Downstream Applications

Total RNA purified with the simplyRNA Tissue Kit or Cells Kit is suitable for use in downstream applications. For more information on downstream applications, see the Promega *Protocols and Applications Guide* (1) and the *RNA Analysis Notebook* (2), available online at **www.promega.com/resources**/

**Note:** For all downstream applications, it is essential that you continue to protect your RNA samples from RNases. Continue to wear clean gloves and use solutions and centrifuge tubes that are RNase-free.

## 8.C. References

- 1. Protocols and Applications Guide, Online Edition (2011) Promega Corporation.
- 2. RNA Analysis Notebook (2002) Promega Corporation.

# 8.D. Related Products

Instrument Accessories	Size	Cat.#
Maxwell® 16 SEV Hardware Kit	1 each	AS1200
Maxwell® 16 LEV Hardware Kit	1 each	AS1250
Maxwell® 16 LEV Cartridge Rack	1 each	AS1251
Maxwell® 16 High Strength LEV Magnetic Rod and Plunger Bar Adaptor	1 each	SP1070
Standard Elution Volume (SEV) Kits	Size	Cat.#
Maxwell® 16 Blood DNA Purification Kit	48 preps	AS1010
Maxwell <sup>®</sup> 16 Cell DNA Purification Kit	48 preps	AS1020
Maxwell® 16 Tissue DNA Purification Kit	48 preps	AS1030
Maxwell® 16 Total RNA Purification Kit	48 preps	AS1050
Maxwell® 16 Mouse Tail DNA Purification Kit	48 preps	AS1120
DNA IQ™ Reference Sample Kit for Maxwell® 16	48 preps	AS1040
Maxwell® 16 Polyhistidine Protein Purification Kit	48 preps	AS1060
LEV Reagent Kits	Size	Cat.#
Maxwell <sup>®</sup> 16 FFPE Tissue LEV DNA Purification Kit	48 preps	AS1130
Maxwell <sup>®</sup> 16 FFPE Plus LEV DNA Purification Kit	48 preps	AS1135
Maxwell® 16 Cell LEV DNA Purification Kit	48 preps	AS1140
Maxwell® 16 Viral Total Nucleic Acid Purification Kit	48 preps	AS1150
Maxwell® 16 Tissue LEV Total RNA Purification Kit	48 preps	AS1220
Maxwell® 16 Cell LEV Total RNA Purification Kit	48 preps	AS1225
Maxwell® 16 LEV Blood DNA Kit	48 preps	AS1290
Maxwell® 16 Buccal Swab LEV DNA Purification Kit	48 preps	AS1295
DNA IQ™ Casework Pro Kit for Maxwell® 16*	48 preps	AS1240
*Not For Medical Diagnostic Use		

\*Not For Medical Diagnostic Use.

Accessory Products	Size	Cat.#	
MagneSphere® Technology Magnetic Separation Stand (twelve-position)	Iagnetic Separation Stand (twelve-position)0.5ml		
Maxwell <sup>®</sup> 16 LEV Magnet	1 each	AS1261	

#### 8.E. Summary of Changes

The following changes were made to the 8/17 revision of this document:

1. Added "to purify RNA from fresh or frozen mammalian cells or fresh or frozen mammalian tissue samples, respectively" to the end of the first sentence of Section 1.

<sup>(a)</sup>U.S. Pat. No. 6,855,499, European Pat. Nos. 1368629, 2090655 and 2363476, Japanese Pat. No. 4399164 and other patents.

© 2011, 2014, 2017 Promega Corporation. All Rights Reserved.

MagneSphere and Maxwell are registered trademarks of Promega Corporation. DNA IQ is a trademark of Promega Corporation.

RNA/ater is a registered trademark of Applied Biosystems. Tissue-Tearor is a trademark of BioSpec Products, Inc.

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.