

# SV Total RNA Isolation System

INSTRUCTIONS FOR USE OF PRODUCTS Z3100, Z3101 AND Z3105.

Quick  
PROTOCOL

## Spin Protocol

1. Place **175µl RNA Lysis Buffer (RLA)** (+ BME) in an autoclaved tube.
2. Prepare sample for lysis.
3. Immediately place sample into **Lysis Buffer**. Mix thoroughly by inversion.  
**Note:** Ensure proper ratio of Lysis Buffer to sample. See Table 1 of the standard protocol.\*
4. Add **350µl RNA Dilution Buffer (RDA, blue)**. Mix by inverting 3–4 times.  
**Note:** Refer to the appropriate lysate preparation section in the Technical Manual #TM048 to determine whether the sample should be heated at 70°C for 3 minutes.
5. Centrifuge for 10 minutes. Transfer the cleared lysate to a fresh tube.
6. Add **200µl 95% ethanol** to cleared lysate and mix well (pipet).

————— **The Spin and Vacuum Protocols are identical up to this point.** —————

7. Transfer mixture to Spin Basket Assembly and centrifuge for 1 minute. Discard eluate.
8. Add **600µl of RNA Wash Solution (RWA)** (+ ethanol). Centrifuge for 1 minute and discard the eluate.
9. Prepare **DNase incubation mix** using the table below:

Solution	Volume	×	Number of Preps	=	Total
Yellow Core Buffer	40µl				
MnCl <sub>2</sub> , 0.09M	5µl				
DNase I	5µl				

*Mix gently (pipet); do **not** vortex.*

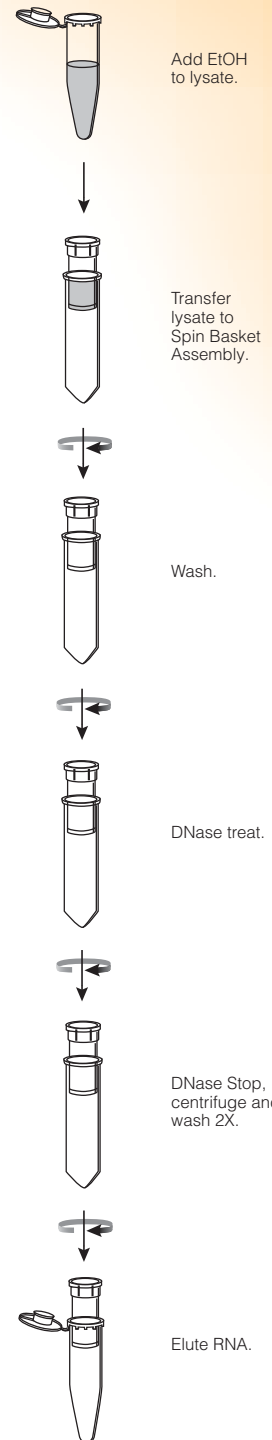
10. Apply **50µl** of DNase mix to membrane. Incubate at RT for 15 minutes.
11. Add **200µl DNase Stop Solution (DSA)** (+ ethanol) and centrifuge for 1 minute.
12. Add **600µl RNA Wash Solution (RWA)**; centrifuge for 1 minute. Empty.
13. Add **250µl RNA Wash Solution (RWA)**; centrifuge for 2 minutes. Transfer Spin Basket to Elution Tube.
14. Add **100µl Nuclease-Free Water** to membrane. Centrifuge for 1 minute to elute the RNA and store at –70°C.

RT: room temperature  
Centrifugation: 12,000–14,000 × *g* (at RT)

\*Additional protocol information is available in Technical Manual #TM048, available online at: [www.promega.com](http://www.promega.com)

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA. Revised 3/09  
Part #9FB025

# SV Total RNA Isolation System

INSTRUCTIONS FOR USE OF PRODUCTS Z3100, Z3101 AND Z3105.

Quick  
PROTOCOL

## Vacuum Protocol

**Note:** For the Vacuum Protocol, follow Steps 1–6 of the Spin Protocol.

- Attach Vacuum Adapter with Luer-Lok® fitting to one manifold port. Gently press SV RNA Spin Basket into adapter and transfer mixture to Spin Basket. Apply vacuum. **Note:** Label Collection Tube and save for Step 13.
- Add **900µl RNA Wash Solution (RWA)**. Apply vacuum until solution has passed through. Stop vacuum source and open unused port to vent manifold. *Release all vacuum pressure before continuing!*
- Prepare **DNase incubation mix** using the table below:

Solution	Volume	×	Number of Preps	=	Total
Yellow Core Buffer	40µl				
MnCl <sub>2</sub> , 0.09M	5µl				
DNase I	5µl				

Mix gently (pipet); do **not** vortex.

- Apply **50µl** of DNase incubation mix to membrane. Incubate at RT for 15 minutes.
- Add **200µl DNase Stop Solution (DSA)** (+ ethanol) to Spin Basket. Close open port and apply vacuum.
- Add **900µl RNA Wash Solution (RWA)**. Repeat wash.
- Release vacuum pressure. Place Spin Basket in Collection Tube (from Step 7). Centrifuge Spin Basket/Collection Tube for 1 minute.
- Transfer Spin Basket to Elution Tube, add **100µl Nuclease-Free Water** and centrifuge for 1 minute. Store purified RNA at  $-70^{\circ}\text{C}$ .

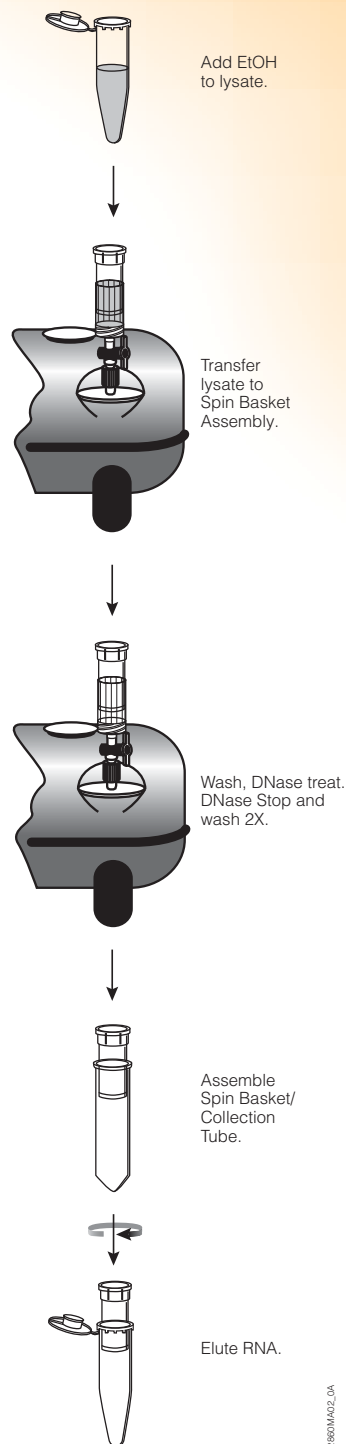
RT: room temperature

Centrifugation:  $12,000\text{--}14,000 \times g$  (at RT)

Additional protocol information is available in Technical Manual #TM048, available online at: [www.promega.com](http://www.promega.com)

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA. Revised 3/09.  
Part #9FB025

© 2000–2009 Promega Corporation. All Rights Reserved.