### QUICK REFERENCE

**Pub. No.** MAN0019349 **Rev.** A.0

Contents Catalog Number 12183020

| Components                              | Amount  | Storage          |  |
|---|---------|------------------|--|
| Lysis Buffer                            | 20 mL   |                  |  |
| Wash Buffer I                           | 10 mL   |                  |  |
| Wash Buffer II                          | 4 mL    |                  |  |
| RNase-free Water                        | 3 mL    | Room temperature |  |
| Spin Cartridges (with collection tubes) | 10 each |                  |  |
| Collection Tubes                        | 10 each |                  |  |
| Recovery Buffer                         | 10 each |                  |  |

### Product description

- The PureLink<sup>™</sup> RNA Mini Kit provides a simple, reliable, and rapid method for isolating high-quality total RNA from a wide variety of samples, including animal and plant cells and tissue, blood, bacteria, yeast, and liquid samples. The purified total RNA is suitable for use in a variety of downstream applications.
- Isolate up to 1 mg of nucleic acid.

## Required materials

2–mercaptoethanol

Online

resources

- 100% ethanol
- (Optional) PureLink<sup>™</sup> DNase Set (Cat. No. 12185010)
- Microcentrifuge capable of 12,000 × *g*
- 1.5 mL RNase-free microcentrifuge tubes
- RNase-free pipet tips

 Visit our product pages for protocols, safety, and additional product information.

- Go online to view related PureLink<sup>™</sup> products.
- For support, visit thermofisher.com/support.

#### Before first use of the kit

#### Prepare Wash Buffer II

Add 16 mL 96–100% ethanol to Wash Buffer II. Check the box on the Wash Buffer II label to indicate that ethanol was added. Store Wash Buffer II with ethanol at room temperature.

#### (Optional) Prepare PureLink<sup>™</sup> DNase Mixture

Add the following components to a clean, RNase-free microcentrifuge tube. Prepare 80 µL for each sample to be processed. Store PureLink<sup>™</sup> DNase Mixture at -20°C.

| Component                   | Volume |
|-----------------------------|--------|
| 10X DNase I Reaction Buffer | 8 µL   |
| Resuspended DNase (~3U/µL)  | 10 µL  |
| RNase free water            | 62 μL  |

### Before each use of the kit

#### Prepare fresh Lysis Buffer

Add 10  $\mu L$  of 2–mercaptoe thanol for each 1 mL of Lysis Buffer needed for the purification procedure in a RN ase-free tube.

Use 1 volume of Lysis Buffer for every 0.2 mL of whole blood (or fraction thereof).

## Troubleshooting

For detailed troubleshooting instructions see the PureLink<sup>™</sup> RNA Mini Kit User Guide at thermofisher.com or contact Technical Support.

# Limited product warranty and licensing information

**Disclaimer:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Manufacturer: Life Technologies Corporation | 2130 Woodward Street | Austin, TX 78744

O2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



## Protocol for purification of RNA from liquid samples

This protocol describes how to purify or desalt one sample of total RNA from  $\leq 1.2$  mL of liquid sample (e.g., enzymatic reaction, upper phase from TRIzol<sup>™</sup> reagent protocol in ethanol) on one column using the PureLink<sup>™</sup> RNA Mini Kit. For detailed instructions see the PureLink<sup>™</sup> RNA Mini Kit User Guide at thermofisher.com or contact Technical Support.

## Important guidelines

• Use proper RNA handling techniques when working with RNA.

| Step |  | <u>כ</u>         | Action  |
|------|--|------------------|---|
| 1    |  | Add Lysis Buffer | <ul> <li>a. Add one volume of liquid sample (≤1.2 mL) to a 1.5 mL RNase-free microcentrifuge tube.</li> <li>b. Add one volume of Lysis Buffer with 2-mercaptoethanol to the sample.</li> <li>c. Add one volume of 100% ethanol to the sample.</li> <li>d. Mix by vortexing or pipetting up and down 5 times.</li> </ul>           |
| 2    |  | Bind RNA         | <ul> <li>a. Transfer the sample (≤700 µL) to a Spin Cartridge (with Collection Tube).</li> <li>b. Centrifuge at 12,000 × g for 15 sec at room temperature. Discard the flow-through, and reinsert the Spin Cartridge into the same Collection Tube.</li> <li>c. Repeat steps a-b until the entire sample is processed.</li> </ul> |

**Thermo Fisher** 

SCIENTIFIC

| Step   |          | Action   |   |  |
|--------|----------|--|---|--|
|        |          | No DNase treatment   | On-column DNase treatment   |  |
| 3<br>3 | Wash RNA | <ul> <li>a. (for cytoplasmic RNA extracts) Add 700 µL Wash<br/>Buffer I to the Spin Cartridge.<br/>Centrifuge at 12,000 × g for 15 sec at room<br/>temperature.<br/>Discard the flow-through and the Collection Tube.<br/>Place the Spin Cartridge into a new Collection<br/>Tube.</li> <li>b. Add 500 µL Wash Buffer II with ethanol to the<br/>Spin Cartridge.</li> <li>c. Centrifuge at 12,000 × g for 15 sec at room<br/>temperature.</li> <li>d. Discard the flow-through and reinsert the Spin<br/>Cartridge in the same Collection Tube.</li> <li>e. Repeat steps b-d one more time.</li> </ul> | <ul> <li>a. Add 350 µL Wash Buffer I to the Spin Cartridge.</li> <li>b. Centrifuge at 12,000 × g for 15 sec at room temperature.</li> <li>c. Discard the flow-through and the Collection Tube. Place the Spin Cartridge into a new Collection Tube.</li> <li>d. Add 80 µL PureLink™ DNase Mixture onto the surface of the Spin Cartridge membrane.</li> <li>e. Incubate at room temperature for 15 min.</li> <li>f. Add 350 µL Wash Buffer I to the Spin Cartridge.</li> <li>g. Centrifuge at ~2,600 × g for 5 min at room temperature.</li> <li>h. Discard the flow-through and the Collection Tube. Place the Spin Cartridge into a new Collection Tube.</li> <li>j. Centrifuge at 12,000 × g for 15 sec at room temperature.</li> <li>k. Discard the flow-through and reinsert the Spin Cartridge.</li> <li>j. Centrifuge at 12,000 × g for 15 sec at room temperature.</li> </ul> |  |
|        |          |  | l. Repeat steps I–k one more time.  |  |

ThermoFisher SCIENTIFIC

# Protocol for purification of RNA from liquid samples

| Step |  | ט                                | Action  |  |
|------|--|----------------------------------|---|--|
| 4    |  | Elute RNA                        | <ul> <li>a. Centrifuge the Spin Cartridge with Collection Tube at 12,000 × g for 1 min at room temperature.</li> <li>b. Discard the Collection Tube and insert the Spin Cartridge into a Recovery Tube.</li> <li>c. Add 30 µL to 3 × 100 µL RNase-Free Water to the center of the Spin Cartridge.</li> <li>d. Incubate at room temperature for 1 min.</li> <li>e. Centrifuge at 12,000 × g for 2 min at room temperature.<br/>Note: Collect all eluates into the same tube when performing serial elution.</li> </ul> |  |
| 5    |  | Analyze RNA yield and<br>quality | Determine the quantity and quality of the purified total RNA using any of the following techniques (See the PureLink™ RNA Mini Kit User Guide for details).<br>• UV absorbance at 260 nm<br>• Fluorescence microplate reader with Quant-iT <sup>™</sup> RiboGreen <sup>™</sup> RNA Assay Kit  |  |
| 6    |  | Store RNA                        | <ul> <li>Keep purified RNA on ice if using the RNA within a few hours of isolation.</li> <li>Store purified RNA at -80°C or long-term storage.</li> </ul>   |  |

24 April 2020