# PureLink<sup>™</sup> miRNA Isolation Kit USER GUIDE

## For rapid, efficient purification of small RNA molecules

Catalog Number K157001 Document Part Number 250753 Publication Number MAN0000474 Revision A.0



The information in this guide is subject to change without notice.

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.

#### Revision history: Pub. No. MAN0000474

Revision	Date	Description
A.0	13 December 2016	Updated to latest branding guidelines.
2.0	27 September 2012	Baseline for revisions.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2016 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

Product information	5
Product description	5
Content and storage	. 5
Required materials not supplied	6
Methods	. 7
Procedural guidelines	7
Before you begin	. 7
Prepare lysates	. 7
Prepare mammalian cell lysates	. 8
Prepare tissue lysates	. 8
Prepare plant lysates	. 8
Prepare yeast lysates	9
Prepare bacterial lysates	9
Prepare lysates from large sample amount	10
Prepare lysates from dicing reaction	10
Isolate RNA	11
Bind the RNA to the membrane	11
Wash the RNA on the membrane	11
Elute the RNA	11
Determine RNA quality and quantity	12
Estimate RNA quantity	12
Analysis of the RNA quality	12
APPENDIX A Troubleshooting	13
APPENDIX B Expected results	14
Results	14
Expected yields	15

APPENDIX C	One-column protocol	16
Bind the RNA to th	ne membrane	16
	the membrane	
Elute the RNA		17
Precipitate the RN	ΙΑ	17
APPENDIX D	Ordering information	18
Additional produc	ts	18
APPENDIX E	Safety	19
Chemical safety .		. 20
Biological hazard	safety	. 21
Documentatio	n and support	22
Customer and tec	hnical support	22
Limited product w	arranty	22
References		23



# **Product information**

**IMPORTANT!** Before using this product, read and understand the information in the "Safety" appendix in this document.

### **Product description**

The Invitrogen<sup>™</sup> PureLink<sup>™</sup> miRNA Isolation Kit allows for isolation of small RNA molecules from biological samples for functional analysis. The kit is specifically designed to isolate high-quality small RNA molecules including tRNA, 5S rRNA, 5.8S rRNA, and regulatory RNA molecules such as microRNA (miRNA) and short interfering RNA (siRNA). PureLink<sup>™</sup> miRNA Isolation Kit uses silica-based membranes to bind the RNA, ensuring high yield and purity and minimal genomic DNA contamination. The isolated RNA using the PureLink<sup>™</sup> miRNA Isolation Kit is suitable for use in Northern blotting and microarray analysis.

This document described RNA isolation from bacteria, yeast, plant, mammalian cells, tissues, and virus.

#### **Content and storage**

Sufficient reagents are included in the kit to perform 25 reactions.

Contents	Amount	Storage
inding Buffer (L3) 45 mL		
Wash Buffer (W5)	10 mL	
Buffer (W4) for single column purification	3 mL	
Sterile, RNase-free Water	5 mL	15°C to 30°C
Spin Cartridges with Collection Tubes	50 cartridges	
Wash Tubes (2 mL)	50 tubes	
Recovery Tubes (1.7 mL)	50 tubes	

Table 1PureLink<sup>™</sup> miRNA Isolation Kit (Cat. No. K157001)



## **Required materials not supplied**

Unless otherwise indicated, all materials are available through **thermofisher.com**. MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

Item	Source		
Equipment			
Benchtop microcentrifuge	MLS		
Mircocentrifuge capable of reaching >10,000 × $g$	MLS		
Laboratory mixer (vortex or equivalent)	MLS		
Homogenizer (for viscous or debris-containing samples only)	12183026		
Homogenizer or tissue grinder (for plant and tissue lysates only)	MLS		
Water bath or heat block at 30°C (for yeast lysate only)			
Tubes, plates, and accessories			
Disposable, individually wrapped, sterile plasticware	MLS		
Sterile, DNase-free microcentrifuge tubes, 1.5 mL	MLS		
Multichannel pipettes	MLS		
Aerosol-resistant pipette tips	MLS		
Reagents			
Ethanol, 70%	MLS		
Ethanol, 96–100%	MLS		
TE Buffer (10 mM Tris-HCl, pH 8.0, 0.1 mM EDTA)	MLS		
Lysozyme (for bacterial lysate only)	MLS		
5% SDS solution (for bacterial lysate only)	MLS		
Zymolase or lyticase (for yeast lysate only)	MLS		
TRIzol <sup>™</sup> Reagent (for large sample amount only)	15596026		
Chloroform (for large sample amount only)	MLS		
UltraPure <sup>™</sup> Glycogen, 20 µg/µL (for one-column protocol only)	10814-010		



# Methods

### **Procedural guidelines**

- Perform all steps at room temperature (20–25°C) unless otherwise noted.
- Use disposable, individually wrapped, sterile plasticware.
- Use only sterile, new pipette tips and microcentrifuge tubes.
- Wear latex gloves while handling reagents and RNA samples to prevent RNase contamination.
- Always use proper microbiological aseptic techniques when working with RNA.
- Use RNase *AWAY*<sup>™</sup> Decontamination Reagent (Cat. No. 10328011) or RNase*Zap*<sup>™</sup> RNase Decontamination Solution (Cat. No. AM9780) to remove RNase contamination from surfaces and from non-disposable equipment like homogenizers.

#### Before you begin

- **For bacterial samples:** Prepare 1% lysozyme in TE Buffer and sterilize by filtration the solution.
- Add 40 mL of 96–100% ethanol to 10 mL of Wash Buffer (W5), then store at room temperature.
- For one-column protocol: Add 1.5 mL of RNase-free Water and 1.5 mL of 96–100% ethanol to 10 mL of Wash Buffer (W4), then store at room temperature.

#### **Prepare lysates**

This section describes protocols to prepare lysates from:

- Mammalian cells (see page 8)
- Mammalian tissues (see page 8)
- Plant tissues (see page 8)
- Yeast (see page 9)
- E. coli (see page 9)
- Large sample amount (see page 10)
- Dicing reaction (see page 10)



Prepare	1. Harvest the cells.
mammalian cell lysates	• Adherent cells (up to 1 × 10 <sup>6</sup> cells): Remove the growth medium from the culture wells.
	• <b>Suspension cells</b> (up to $1 \times 10^6$ cells): Centrifuge the culture plate at $250 \times g$ for 5 minutes, then remove the growth medium.
	<b>2.</b> Add 300 μL of Binding Buffer (L3).
	<b>3.</b> Mix well by vortexing or pipetting up and down.
	<b>4.</b> Add 300 μL of 70% ethanol.
	<b>5.</b> Mix well by vortexing.
	Proceed immediately to "Isolate RNA" on page 11.
Prepare tissue	1. Place ~5 mg of minced mammalian tissue in a sterile microcentrifuge tube.
lysates	<ol> <li>Add 300 μL of Binding Buffer (L3).</li> <li>Ensure that the tissue is completely immersed in the buffer.</li> </ol>
	<b>3.</b> Homogenize the tissue using a tissue homogenizer.
	<b>4.</b> Centrifuge the lysate for 5 minutes at $12,000 \times g$ .
	<b>5.</b> Transfer the supernatant to a fresh tube.
	<b>Note:</b> If the lysate is viscous or contains cell debris, clarify the lysate using the Homogenizer (Cat. No. 12183026).
	<b>6.</b> Add 300 μL of 70% ethanol.
	<b>7.</b> Mix well by vortexing.
	Proceed immediately to "Isolate RNA" on page 11.
Prepare plant	1. Prepare plant tissue:
lysates	<ul> <li>Hard tissue: Freeze the plant tissue in liquid nitrogen and grind to a powder.</li> <li>Soft tissue: Cut soft, non-fibrous plant tissue into small pieces.</li> </ul>
	<b>2.</b> Add 300 $\mu$ L of Binding Buffer (L3).
	<b>3.</b> Prepare lysate.
	• <b>Hard tissue:</b> Vortex the ground tissue until the powder is completely resuspended.
	• <b>Soft tissue:</b> Homogenize with a homogenizer or tissue grinder.
	<b>4.</b> Centrifuge the lysate at $12,000 \times g$ for 2 minutes to remove insoluble materials.



	<b>5.</b> Transfer the supernatant to a new microcentrifuge tube.
	<b>Note:</b> If the lysate is viscous or contains cell debris, clarify the lysate using the Homogenizer (Cat. No. 12183026).
	<b>6.</b> Add 300 μL of 70% ethanol.
	<b>7.</b> Mix well by vortexing.
	Proceed immediately to "Isolate RNA" on page 11.
Prepare yeast	If you are using a frozen cell pellet, proceed directly to step 2.
lysates	<b>1.</b> Harvest 1 mL of fresh, log-phase yeast cells (OD <sub>660</sub> =1.04) by centrifugation.
	<b>2.</b> Resuspend the pellet in 70 $\mu$ L of cold TE Buffer.
	<b>3.</b> Add 30 units of zymolase (or lyticase).
	<b>4.</b> Centrifuge the lysate at $12,000 \times g$ for 2 minutes to remove insoluble materials.
	<b>5.</b> Resuspend the pellet in 300 $\mu$ L of Binding Buffer (L3).
	<b>6.</b> Add 210 μL of 96-100 ethanol.
	<b>7.</b> Mix well by vortexing.
	Proceed immediately to "Isolate RNA" on page 11.
Prepare bacterial	If you are using a frozen cell pellet, proceed directly to step 2.
lysates	<b>1.</b> Harvest up to $5 \ge 10^6 E$ . <i>coli</i> cells by centrifugation.
	<b>2.</b> Resuspend the pellet in 90 $\mu$ L of 1% lysozyme in TE Buffer.
	<b>3.</b> Add 1 $\mu$ L of 5% SDS to the lysate and mix thoroughly.
	<b>4.</b> Incubate for 5 minutes.
	<b>5.</b> Add 300 $\mu$ L of Binding Buffer (L3).
	<b>6.</b> Mix well by vortexing.
	<b>7.</b> Add 210 μL of 96-100 ethanol.
	<b>8.</b> Mix well by vortexing.
	Proceed immediately to "Isolate RNA" on page 11.



#### Prepare lysates from large sample amount

**1.** Lyse the samples in TRIzol<sup>™</sup> Reagent.

Sample type	Lysis procedure
Tissues	Homogenize tissue samples in 1 mL of TRIzol <sup>™</sup> Reagent per 50– 100 mg of tissue using a tissue homogenizer.
Adherent cells	<ol> <li>Add 1 mL of TRIzol<sup>™</sup> Reagent per 10 cm<sup>2</sup> directly to the culture dish.</li> <li>Lyse the cells by passing the cell lysate several times through a pipette.</li> </ol>
Suspension Cells	<ol> <li>Harvest cells</li> <li>Harvest and pellet the cells by centrifugation.</li> <li>Add 1 mL of TRIzol<sup>™</sup> Reagent per 5–10 × 10<sup>6</sup> animal, plant, or yeast cells, or per 1 × 10<sup>7</sup> bacterial cells.</li> <li>Lyse the cells by passing the cell lysate several times through a pipette.</li> </ol>

- **2.** Incubate for 5 minutes.
- Add 0.2 mL of chloroform per 1 mL of TRIzol<sup>™</sup> Reagent, then shake the tube vigorously by hands for 15 seconds.
   Do not vortex the tube.

**4.** Centrifuge the sample at  $12,000 \times g$  for 15 minutes at 4°C. The mixture separates into a lower red phenol-chloroform phase, an interphase, and a colorless upper aqueous phase. The volume of the aqueous phase is ~600 µL.

- 5. Transfer ~400  $\mu$ L of the colorless, upper phase containing RNA to a fresh tube.
- **6.** Add 215 μL of 96–100% ethanol.
- **7.** Mix well by vortexing.

Proceed immediately to "Isolate RNA" on page 11.

Prepare lysates from dicing reaction

1. Add 300  $\mu$ L of Binding Buffer (L3) and 300  $\mu$ L of 96–100% ethanol to 300  $\mu$ L of the dicing reaction.

Follow this procedure to prepare samples produced in a dicing reaction.

2. Mix well by vortexing.

Proceed immediately to "Isolate RNA" on page 11.



### **Isolate RNA**

Bind the RNA to the membrane	<b>1.</b> Transfer sample to a Spin Cartridge with Collection Tube.
	<b>2.</b> Centrifuge for 1 minute at $12,000 \times g$ .
	<b>3.</b> Add 700 μL of 96–100% ethanol.
	<b>4.</b> Mix well by vortexing.
	<b>5.</b> Transfer 700 $\mu$ L of the sample to a new Spin Cartridge.
	<b>6.</b> Centrifuge for 1 minute at $12,000 \times g$ .
	<b>7.</b> Transfer remaining sample to the Spin Cartridge from step 5.
	<b>8.</b> Discard the flow through, then place the Spin Cartridge into the Collection Tube.
Wash the RNA on	1. Add 500 $\mu$ L of Wash Buffer to the Spin Cartridge.
the membrane	<b>2.</b> Centrifuge for 1 minute at $12,000 \times g$ .
	<b>3.</b> Repeat step 1-step 2.
	<b>4.</b> Discard the flow through, then place the Spin Cartridge into a Wash Tube.
	<b>5.</b> Centrifuge for 2–3 minutes at maximum speed.
	<b>6.</b> Discard the Wash Tube.
Elute the RNA	1. Place the Spin Cartridge into a clean 1.7–mL Recovery Tube.
	<b>2.</b> Add 50–100 $\mu$ L of Sterile, RNase-free Water to the center of the cartridge.
	<b>3.</b> Incubate for 1 minute.
	<b>4.</b> Centrifuge for 1 minute at maximum speed.
	<b>5.</b> Remove and discard the Spin Cartridge. The Recovery Tube contains the purified RNA.

Proceed immediately to downstream application or store the purified RNA at  $-80^{\circ}$ C for further use.

# Determine RNA quality and quantity

Estimate RNA quantity	Alternatively, the purified RNA can be quantified using Quant-iT <sup>™</sup> RNA Assay Kit (Cat. No. Q33140).		
	<b>1.</b> Dilute an aliquot of the total RNA sample in 10 mM Tris-HCl, pH 7.0.		
	<b>2.</b> Mix well, then transferred to a cuvette (1-cm path length).		
	<b>Note:</b> The RNA must be in a neutral buffer to accurately measure the UV absorbance.		
	<b>3</b> . Determine the OD <sub>260</sub> of the solution using a spectrophotometer blanked against 10 mM Tris-HCl, pH 7.0.		
	4. Calculate the amount of total RNA using the following formula:		
	Total RNA ( $\mu$ g) = OD <sub>260</sub> × [40 $\mu$ g/(1 OD <sub>260</sub> × 1 mL)] × dilution factor × total sample volume (mL)		
Analysis of the RNA quality	Typically, RNA isolated using the PureLink <sup>TM</sup> miRNA Isolation Kit has an $OD_{260/280}$ of >1.8 when samples are diluted in Tris-HCl (pH 7.5). An $OD_{260/A280}$ of >1.8 indicates that RNA is reasonably clean of proteins and other UV chromophores (heme, chlorophyll, etc.) that could either interfere with downstream applications or negatively affect the stability of the stored RNA.		
	To visualize small RNA molecules, you can perform a denaturing gel electrophoresis using 10% or 15% TBE acrylamide gels (see page 14 for an example).		



# Troubleshooting

Observation	Possible cause	Recommended action
The RNA yield is low	The lysis is incomplete or too much sample has clogged the filter.	Decrease the amount of starting material used.
		For tissues, cut the tissue into smaller pieces and ensure that the tissue is completely immersed in Binding Buffer (L3) to obtain optimal lysis.
		If the cartridge is clogged, decrease the sample volume used or load the sample on 2 Spin Cartridges.
	The sample has a low RNA content.	Various tissues have different small RNA content and some tissue may not contain any small RNA at detectable levels.
	The flow through from the first Spin Cartridge was discarded.	Do not discard the flow through from the first cartridge. The flow through contains small RNA molecules.
	The binding conditions were incorrect.	For efficient binding of small RNA molecules, always add ethanol to the flow through to a final concentration of 70%.
	Ethanol was not added to Wash Buffer (W5).	Be sure to add 96–100% ethanol to Wash Buffer (W5) as described in "Before you begin" on page 7.
	The elution conditions were incorrect.	Add Sterile, RNase-free Water to the center of the Spin Cartridge, then incubate for 1 minute before centrifugation.
	The RNA was quantified in water.	Be sure the RNA quantitation using UV absorbance is performed with 10 mM Tris-HCl, pH 7.0 to accurately measure the UV absorbance.
The RNA is degraded	The RNA is contaminated by RNase.	Follow "Procedural guidelines" on page 7.
	The starting materials are of poor quality.	Always use fresh samples or samples frozen at –80°C. For lysis, process the sample quickly to avoid degradation.
The RNA is contaminated by genomic DNA	A large sample amount was loaded onto the first Spin	Follow "Prepare lysates from large sample amount" on page 10.
	Cartridge.	Perform DNase I digestion on the RNA sample after elution to remove genomic DNA contamination.
The RNA is contaminated by total RNA	A large sample amount was loaded onto the first Spin Cartridge.	Follow "Prepare lysates from large sample amount" on page 10.



# **Expected results**

### Results

An example of small RNA molecules isolated from various samples is shown in the following figure.



Figure 1 Small RNA molecules from various samples isolated using PureLink<sup>™</sup> miRNA Isolation Kit

Samples (10  $\mu$ L eluate) were analyzed on a Novex TBE-Urea 15% Gel (Cat. No. EC6885BOX) and RNA bands were visualized with ethidium bromide staining after electrophoresis.

- (1) 1 µL 10 bp DNA Ladder (5) Mouse liver
- 2 HeLa cells
- 3 293F cells
- ④ Rat spleen

- 6 Spinach
  - ⑦ MaV yeast cells
  - (8) *E. coli* cells

### **Expected yields**

The yield of small RNA molecules obtained from various samples using PureLink<sup>TM</sup> miRNA Isolation Kit is listed in the following table. The RNA quantitation was performed with the Quant-iT<sup>TM</sup> RNA Assay Kit (Cat. No. Q33140).

Table 2	Yield of total RNA obtained from various sources	

Source	Amount	small RNA yield
HeLa cells	1 × 10 <sup>6</sup> cells	1.29 µg
293F cells	1 × 10 <sup>6</sup> cells	1.95 µg
Mouse liver	5 mg	0.71 µg
Rat spleen	5 mg	1.39 µg
Spinach	60 mg	1.64 µg
Yeast ( <i>S. cerevisae</i> )	1 × 10 <sup>6</sup> cells	6.21 µg
Bacteria ( <i>E. coli</i> )	$2 \times 10^{6}$ cells	0.55 µg



# One-column protocol

The one-column purification protocol is only applicable to samples obtained from dicing reaction and is designed for use with up to 8  $\mu$ g of RNA in 50- $\mu$ L dicing reaction as starting material.

- To process <8 μg in <50-μL dicing reaction, increase the volume of the starting material to 50 μL.
- To process >8 µg in >50-µL dicing reaction, adjust the buffer volumes accordingly and perform multiple loadings of the sample onto the Spin Cartridge.

The one-column protocol is not recommended with samples prepared from tissues or cells as the amount of sample used for purification may clog the Spin Cartridges, and the volume and amount of buffers used for wash and elution steps are not sufficient to produce high-quality RNA.

### Bind the RNA to the membrane

- 1. Add 150 µL of Binding Buffer (L3) to the 50-µL dicing reaction, them mix well.
- **2.** Add 600  $\mu$ L of 96–100% ethanol, then mix well.
- **3.** Transfer 700  $\mu$ L of the sample to a Spin Cartridge in a Collection Tube.
- **4.** Centrifuge for 1 minute at  $12,000 \times g$ , then discard the flow through.

#### Wash the RNA on the membrane

- 1. Add 500 µL of Wash Buffer (W5) to the Spin Cartridge.
- **2.** Centrifuge for 1 minute at  $12,000 \times g$ .
- 3. Repeat step 1-step 2.
- 4. Discard the Collection Tube, then place the Spin Cartridge in a Wash Tube.
- **5.** Centrifuge for 1 minute at  $12,000 \times g$ , then discard the Wash Tube.

### **Elute the RNA**

- 1. Place the Spin Cartridge in a Recovery Tube, then add 100  $\mu$ L of Buffer (W4) for single-column purification.
- 2. Incubate for 1 minute.
- **3.** Centrifuge for 1 minute at  $12,000 \times g$ , then discard the Spin Cartridge. The eluate contains the RNA.

### **Precipitate the RNA**

- 1. Add 200  $\mu L$  of cold 96–100% ethanol and 1  $\mu L$  of glycogen solution (20  $\mu g/\mu L)$  to the eluate.
- **2.** Mix well and incubate for 15 minutes at –20°C.
- 3. Centrifuge for 15 minutes at maximum speed at 4°C.
- 4. Carefully discard the supernatant without disturbing the pellet.
- **5.** Resuspend the pellet in 0.5 mL of cold 70% ethanol.
- 6. Centrifuge for 10 minutes at maximum speed at 4°C.
- 7. Carefully discard the supernatant without disturbing the pellet.
- **8.** Air-dry the pellet for ~5 minutes.
- 9. Resuspend the pellet in 50 µL of Sterile, RNase-free Water.

Proceed immediately to downstream application or store the purified RNA at -80°C for further use.



# Ordering information

# Additional products

The following products	are also available through <b>thermofisher.com</b> .
------------------------	--

Item	Quantity	Cat. No.
RNase $AW\!AY^{M}$ Decontamination Reagent	250 mL	10328011
RNase <i>Zap</i> <sup>™</sup> RNase Decontamination Solution	250 mL	AM9780
UltraPure <sup>™</sup> DEPC-treated Water	1 L	750023
UltraPure <sup>™</sup> DNase/RNase-Free Distilled Water	500 mL	10977-015
Quant-iT <sup>™</sup> RNA Assay Kit	1 kit	Q33140
Novex TBE-Urea Gels, 15% (1.0 mm, 10 well)	1 box	EC6885BOX
Novex TBE-Urea Gels, 10% (1.0 mm, 10 well)	1 box	EC6875BOX
Novex TBE-Urea Sample Buffer (2X)	10 mL	LC6876
Novex TBE Running Buffer (5X)	1 L	LC6675
Homogenizer	50 each	12183026
10 bp DNA Ladder	50 µg	10821-015
PureLink <sup>™</sup> RNA Mini Kit	50 reactions	12183018A
TrackIt <sup>™</sup> 10 bp DNA Ladder	20 applications	10488-019
UltraPure <sup>™</sup> Glycogen	100 µL	10814-010
TRIzol <sup>™</sup> Reagent	100 mL	15596026

# Safety





- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see the "Documentation and Support" section in this document.

### **Chemical safety**



**WARNING!** GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

### **Biological hazard safety**



**WARNING!** BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:
- www.cdc.gov/biosafety/publications/bmbl5/BMBL.pdf
  World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:
  www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf

# **Documentation and support**

### **Customer and technical support**

Visit **thermofisher.com/support** for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

#### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **www.thermofisher.com/us/en/home/global/ terms-and-conditions.html**. If you have any questions, please contact Life Technologies at **www.thermofisher.com/support**.

# References

Elbashir, S. M., Harborth, J., Weber, K., and Tuschl, T. 2002. *Analysis of Gene Function in Somatic Mammalian Cells Using Small Interfering RNAs*. Methods 26, 199-213.

Lim, L. P., Glasner, M. E., Yekta, S., Burge, C. B., Bartel, D. P. 2003. *Vertebrate microRNA Genes*. Science 299, 1540.

Liu, C. G., Calin, G. A., Meloon, B., Gamliel, N., Sevignani, C., Ferracin, M., Dumitru, C. D., Shimizu, M., Zupo, S., Dono, M., Alder, H., Bullrich, F., Negrini, M., and Croce, C. M. 2004. *An Oligonucleotide Microchip for Genome-wide microRNA Profiling in Human and Mouse Tissues*. Proc. Natl. Acad. Sci. USA 101, 9740-9744.



