TRIzol[™] Plus RNA Purification Kit

Catalog Numbers 12183555

Doc. Part No. 25-0915 Pub. No. MAN0000561 Rev. A.0

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product information

Invitrogen[™] TRIzol[™] Plus RNA Purification 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, bacteria, and yeast. The kit utilizes the strong lysis capability of TRIzol[™] Reagent, followed by a convenient and time-saving silicacartridge purification protocol from the PureLink[™] RNA Mini Kit, to purify ultrapure total RNA within an hour, even from difficult samples such as fibrous tissue.

TRIzol[™] Reagent is a monophasic solution of phenol, guanidine isothiocyanate, and other proprietary components which facilitate the isolation of a variety of RNA species of large or small molecular size. TRIzol[™] Reagent maintains the integrity of the RNA due to highly effective inhibition of RNase activity while disrupting cells and dissolving cell components during sample homogenization or lysis.

TRIzol[™] Plus RNA Purification Kit can also be used with Phasemaker[™] Tubes to isolate RNA. Refer to *TRIzol[™] Plus RNA Purification Kit and Phasemaker[™] Tubes Complete System User Guide* (MAN0016165) for the full protocol.

Contents and storage

Reagents provided in the kit are sufficient for 50 RNA isolations.

Table 1 TRIzol[™] Plus RNA Purification Kit (Cat. No. 12183555)

Contents	Amount	Storage
TRIzol [™] Reagent	100 mL	15-30°C
PureLink [™] RNA Mini Kit ^[1]	1 kit	15-50 C

^[1] See Table 2 for kit contents.

Table 2 PureLink[™] RNA Mini Kit (Cat. No. 12183018A)

Contents	Amount	Storage
Spin cartridges (with collection tubes)	50 cartridges	
Collection tubes	50 tubes	
Wash Buffer I	50 mL	45,0000
Wash Buffer II	15 mL	15–30°C
RNase-free water	15.5 mL	
Recovery tubes	50 tubes	
Lysis Buffer ^[1]	125 mL	

^[1] Not used in this protocol.

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			
Homogenizer, or equivalent	12183026		
Tubes and tips			
1.5-mL RNase-free microcentrifuge tubes	MLS		
RNase-free pipette tips	MLS		
Reagents			
Chloroform, or 4-bromoanisole	MLS		
Ethanol, molecular grade	MLS		

Input sample requirements

IMPORTANT! Perform RNA isolation immediately after sample collection or quick-freeze samples immediately after collection and store at -80°C or in liquid nitrogen until RNA isolation.

Sample type	Starting material per 1 mL of TRIzol™ Reagent
Tissues ^[1]	50–100 mg of tissue
Cells grown in monolayer	1 × 10 ⁵ –1 × 10 ⁷ cells grown in monolayer in a 3.5–cm culture dish (10 cm ²)
Cells grown in suspension	$5-10 \times 10^6$ cells from animal, plant, or yeasty origin or 1×10^7 cells of bacterial origin

 Fresh tissues or tissues stored in RNA*later*[™] Stabilization Solution (Cat. No. AM7020).

Procedural guidelines

- Perform all steps at room temperature (20–25°C) unless otherwise noted.
- Use disposable, individually wrapped, sterile plastic ware and sterile, disposable RNase-free pipettes, pipette tips, and tubes.
- Wear disposable gloves while handling reagents and RNA samples to prevent RNase contamination from the surface of the skin; change gloves frequently, particularly as the protocol progresses from crude extracts to more purified materials.
- Always use proper microbiological aseptic techniques when working with RNA.
- Use RNaseZap[™] RNase Decontamination Solution (Cat. no. AM9780) to remove RNase contamination from work surfaces and non-disposable items such as centrifuges and pipettes used during purification.
- Always maintain a ratio of 10:1 between the volume of TRIzol[™] Plus RNA Purification Kit and the mass of the sample.
- The maximum RNA binding capacity of the spin cartridge is ~1 mg. If you are processing samples that contain more than 1 mg of total RNA, divide the sample into aliquots containing <1 mg total RNA for each spin cartridge used.



• Use clear polypropylene disposable tubes when working with <2 mL volumes of TRIzol[™] Reagent. For larger volumes, use glass or polypropylene tubes, and ensure that the tubes can withstand centrifugation at 12,000 × *g* with TRIzol[™] Reagent and chloroform. Do not use tubes that leak or crack.

Before first use of the kit

- 1. Add 60 mL of 96–100% ethanol to Wash Buffer II, then mix thoroughly.
- 2. Check the box on the Wash Buffer II label to indicate that ethanol was added, then store at room temperature.

Isolate RNA

1	Lyse samples and separate phases	a.	Lyse and homogenize samples in TRIzol [™] Reagent according to your starting material. • Tissues:
			Add 1 mL of TRIzol [™] Reagent per 50–100 mg of tissue to the sample and homogenize using a homogenizer.
			Cell grown in monolayer:
			a. Remove growth media.
			b. Add 1 mL of TRIzol ^{T} Reagent per 1 × 10 ⁵ -10 ⁷ cells directly to the culture dish to lyse the cells.
			c. Pipet the lysate up and down several times to homogenize.
			Cells grown in suspension:
			 a. Pellet the cells by centrifugation and discard the supernatant. b. Add 0.75 mL of TRIzol[™] Reagent per 0.25 mL of sample (5–10 × 10⁶ cells from animal, plant, or yeasty origin or 1 ×10⁷ cells of bacterial origin) to the pellet.
			Note: Do not wash cells before addition of TRIzol [™] Reagent to avoid mRNA degradation. c. Pipet the lysate up and down several times to homogenize.
			STOPPING POINT Samples can be stored at 4° C overnight or at -20° C for up to a year.
		b.	(<i>Optional</i>) If samples have a high fat content, centrifuge the lysate for 5 minutes at $12,000 \times g$ at $4-10^{\circ}$ C, then transfer the clear supernatant to a new tube.
		c.	Incubate for 5 minutes to permit complete dissociation of the nucleoproteins complex.
		d.	Add 0.2 mL of chloroform or 50 µL of 4-bromoanisole per 1 mL of TRIzol [™] Reagent used for lysis, then securely cap the tube.
		e.	Incubate for 2–3 minutes.
		f.	Centrifuge the sample for 15 minutes at 12,000 × g at 4°C.
			The mixture separates into a lower red phenol-chloroform, and interphase, and a colorless upper aqueous phase.
		g.	Transfer ~600 μ L of the colorless, upper aqueous phase containing the RNA to a new tube.
		h.	Add an equal volume of 70% ethanol, then mix well by vortexing.
		i.	Invert the tube to disperse any visible precipitate that may form after adding ethanol.
2	Bind the RNA to the	a.	Transfer up to 700 μ L of the sample to a spin cartridge (with collection tube).
4	membrane	b.	Centrifuge at $12,000 \times g$ for 15 seconds.
		c.	Discard the flow-through, then reinsert the spin cartridge into the same collection tube.
		d.	Repeat step 2a-step 2c until the entire sample has been processed.
		e.	(<i>Optional</i>) If your downstream application required DNA-free total RNA, perform on-column DNase treatment (see <i>PureLink[™] RNA Mini Kit User Guide</i> (Pub. No. MAN0000406).
3	Wash the RNA on the	a.	Add 700 μ L of Wash Buffer I to the spin cartridge.
9	membrane	b.	Centrifuge at $12,000 \times g$ for 15 seconds.
		с.	Discard the flow-through, then reinsert the spin cartridge into the same collection tube.
			Add 500 μ L of Wash Buffer II to the spin cartridge.
			Centrifuge at 12,000 × g for 15 seconds.
			Discard the flow-through, then reinsert the spin cartridge into the same collection tube.
		g.	Repeat step 3d-step 3f once.
4	Elute the RNA	a.	Centrifuge at $12,000 \times g$ for 1 minute to dry the membrane.
-		b.	Discard the collection tube, then insert the spin cartridge into a recovery tube.
		c.	Add 30 μ L-3 × 100 μ L (3 sequential elutions with 100 μ L each) of RNase-free water to the center of the spin cartridge.
			Note: If you are performing sequential elutions, collect all eluates in the same tube.
			Incubate for 1 minute.
			Centrifuge at >12,000 × g for 2 minutes.
			Discard the spin cartridge.
			e recovery tube contains the purified total RNA.
			Store the purified RNA on ice if used within a few hours. For long-term storage, store the purified RNA at -80°C.
			If highly pure RNA without genomic DNA contamination is required, perform DNase I treatment after purification (see <i>PureLink</i> [™] <i>RNA Mini Kit User Guide</i> (Pub. No. MAN0000406).

Elute the RNA (continued) • You can determine the quality and quantity of the purified RNA using UV absorbance at 260 nm or with the Quant-iT[™] RNA Assay Kit (Cat. No. Q33140).

Troubleshooting

Observation	Possible cause	Recommended action
A lower RNA yield is observed	The samples were incompletely homogenized or lysed.	Use the appropriate method for lysate preparation based on your starting material (see "Input sample requirements" on page 1).
		Decrease the amount of starting material.
		Cut tissue samples into smaller pieces and ensure the tissue is completely immersed in TRIzol [™] Reagent to achieve total lysis.
	The starting material is of poor quality.	Use fresh samples immediately after collection or freeze sample at – 80°C or in liquid nitrogen until further use.
	The spin cartridge is clogged.	Clear lysate and remove any particulate or viscous material by centrifugation, and use only the supernatant for subsequent loading onto the spin cartridge.
	Ethanol was not added to Wash Buffer II .	See "Before first use of the kit" on page 2.
	The elution conditions are incorrect.	Add RNase-free water (30 $\mu\text{L}\mathcal{L}\mathcal{-}3$ × 100 μL) and incubate for 1 minute before centrifugation.
		To recover more RNA, use up to 3 sequential elutions of 100 μL each.
The RNA is degraded	The RNA is contaminated with RNase.	Follow "Procedural guidelines" on page 1 to prevent RNase contamination.
	The samples were improperly handled from collection until lysis.	Perform RNA isolation immediately after sample collection or quick- freeze samples immediately after collection and store at –80°C or in liquid nitrogen until RNA isolation.
Downstream reactions are inhibited	Ethanol is present in the purified RNA.	The spin cartridge membrane has to be dry before eluting the RNA, as traces of ethanol from Wash Buffer II can inhibit downstream enzymatic reactions.
		1. Discard Wash Buffer II flow-through.
		 Insert the spin cartridge into the collection tube. Centrifuge for 1–2 minutes at 12,000 × g to completely dry the membrane.
		4. Proceed with RNA elution.
	Salts are present in the purified RNA.	Always wash with Wash Buffer I before washing with Wash Buffer II.
The RNA $A_{260/280}$ ratio is low	The sample was diluted in water.	Use 10 mM Tris-HCl (pH 7.5) to dilute the samples for absorbance measurements.

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.

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. MAN0000561

Revision	Date	Description
A.0	09 November 2016	Added references to Phasemaker [™] Tubes .
-	29 November 2010	Baseline for revision.

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

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



thermofisher.com