



RNA isolation

User manual

NucleoZOL

This product distributed by Clontech Laboratories, Inc.
A Takara Bio Company
1290 Terra Bella Ave., Mountain View, CA 94043
www.clontech.com

For ordering information
1.800.662.2566
orders@clontech.com

For technical support
1.800.662.2566
tech@clontech.com



For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Clontech Laboratories, Inc. ©2015

July 2015 / Rev. 01

MACHEREY-NAGEL

www.mn-net.com



Table of contents

1	Components	4
1.1	Kit contents	4
1.2	Reagents, consumables, and equipment to be supplied by the user	4
1.3	RNase-free work environment	5
1.4	About this user manual	5
2	Product description	6
2.1	The basic principle	6
2.2	Product specifications	7
2.3	Handling, preparation, and storage of starting materials	8
2.4	RNA reconstitution	8
3	Storage conditions and preparation of working solutions	9
4	Safety instructions	10
5	NucleoZOL protocols	12
5.1	Isolation of small and large RNA in two separate fractions	12
5.2	Isolation of total RNA	15
5.3	Isolation of total RNA in combination with NucleoSpin® RNA kits	18
6	Appendix	20
6.1	Digestion of residual DNA solution	20
6.2	Troubleshooting	21
6.3	Ordering information	22
6.4	Product use restriction / warranty	23

1 Components

1.1 Kit contents

NucleoZOL	
REF	740404.200
NucleoZOL reagent	200 mL

1.2 Reagents, consumables, and equipment to be supplied by the user

Reagents

- RNase-free water
- 75 % ethanol
- 70 % isopropanol
- 100 % isopropanol

Consumables

- 1.5 mL, 2.0 mL or 15 mL centrifuge tubes (depending on the amount of sample to be processed per preparation)
- Sterile RNase-free tips

Equipment

- Manual pipettors
- Vortex mixer
- Centrifuge for microcentrifuge tubes
- Equipment for sample disruption and homogenization
- Personal protection equipment (e.g., lab coat, gloves, goggles)
- Well ventilated working environment
- RNase-free working environment

1.3 RNase-free work environment

The reagent has been tested for functionality. However, an RNase-free working environment is also a critical factor for performing successful RNA isolation and handling. Therefore, general recommendations to avoid RNase contamination should be followed:

- Maintain a separate area, dedicated pipettors, and materials when working with RNA.
- Wear gloves when handling RNA and reagents to avoid contact with skin, which is a source of RNases. Change gloves frequently.
- Use sterile RNase-free plastic tubes. Tubes for lysate preparation and RNA precipitation have to be supplied by the user.
- Keep all kit components sealed when not in use and all tubes tightly closed when possible.

1.4 About this user manual

Please read the detailed protocol if using NucleoZOL for the first time. Experienced users may refer to the short instruction manual.

All technical literature is available on the Internet at www.mn-net.com.

Please contact Technical Service regarding information about changes of to the current user manual compared to the previous revisions.

2 Product description

2.1 The basic principle

NucleoZOL is designed for the isolation of total RNA (small and large RNA) in a single fraction or in separate fractions from a variety of sample materials, such as cells, tissue, and liquids from human or animal origin, plants, yeast, bacteria, viral materials, and other sources.

One of the most important factors during the isolation of RNA is to prevent degradation. First, cells and tissues are lysed and homogenized in NucleoZOL reagent based on guanidinium thiocyanate and phenol. Contaminating molecules such as DNA, polysaccharides, and proteins are precipitated by the addition of water and removed by centrifugation. The NucleoZOL procedure allows the separate isolation of small and large RNA by adding ethanol and isopropanol, respectively. RNA can be reconstituted by RNase-free water. A chloroform-induced phase separation is not necessary for high-quality RNA isolation.

The RNA is ready for use in qRT-PCR, microarrays, RNase protection assays, poly A⁺ isolation, blotting, and other applications.

2.2 Product specifications

Table 1: Product specifications at a glance

Technology	One-phase extraction
Sample material (per mL NucleoZOL)	< 1 x 10 ⁷ cultured cells, bacteria, and yeast, < 100 mg human/animal/plant tissue, < 0.4 mL (viral) fluids
Fragment size	Small RNA (10–200 nt), large RNA (> 200 nt)
Typical yield (total RNA)	Liver: 6–8 µg/mg tissue Kidney, spleen: 3–4 µg/mg tissue Muscle, brain, lung: 0.5–1.5 µg/mg Cultured cells: 4–10 µg/10 ⁶ cells
Typical yield (large RNA)	Liver: 5–7 µg/mg tissue Kidney, spleen: 3–4 µg/mg tissue Muscle, brain, lung: 0.5–1.5 µg/mg Cultured cells: 3–8 µg/10 ⁶ cells
A _{260/280} (total RNA)	1.8–2.1
Typical RIN (RNA integrity number)	> 9
Elution volume	flexible

2.3 Handling, preparation, and storage of starting materials

Sample harvest and RNase inhibition

RNA is not protected against digestion until the sample material is flash frozen or disrupted in the presence of RNase inhibiting or denaturing agents.

Sample harvest methods:

- Use freshly harvested sample for immediate lysis and RNA purification.
- Samples can be stored in NucleoZOL after disruption at -20 °C to -70 °C for up to one year, at 4 °C for up to 24 hours, or up to several hours at room temperature. Frozen samples in NucleoZOL should be thawed slowly before starting with the isolation of RNA.
- Flash freeze sample in liquid N₂ immediately upon harvest and store at -70 °C. Frozen samples are stable up to 6 months. Mortar and pestle can be used to pulverize the sample in a frozen state. Make sure that the sample does not thaw prior to contact with the reagent.
- Samples can be submerged and stored in RNA stabilizing reagents such as RNA/later[®]. Remove excess RNA/later[®] solution from the tissue before processing the sample.

2.4 RNA reconstitution

The precipitated RNA can be dissolved in variable volumes of RNase-free water (see ordering information, section 6.3) to approach an RNA concentration of approximately 1–2 µg/µL for the large RNA fraction and approximately 0.1 µg/µL for the small RNA fraction.

3 Storage conditions and preparation of working solutions

Attention:

NucleoZOL contains phenol (corrosive liquid/poison) and guanidium thiocyanate (irritant). Wear gloves, eye protection, and goggles!

CAUTION: *Read the warning note on the container and MSDS. NucleoZOL contains phenol and guanidinium thiocyanate which CAUSES BURNS and can be fatal. When working with NucleoZOL, use gloves and eye protection (face shield, safety goggles). Do not get the reagent on skin or clothing. Avoid breathing fumes. In case of contact: Immediately flush eyes or skin with a large amount of water for at least 15 minutes and if necessary seek medical attention.*

NucleoZOL can be stored at room temperature (18–25 °C) and is stable for at least one year.

4 Safety instructions


NucleoZOL contains hazardous contents.

Wear gloves and goggles and follow the safety instructions given in this section.

GHS classification

Only harmful features do not need to be labeled with H and P phrases up to 125 mL or 125 g.

Mindergefährliche Eigenschaften müssen bis 125 mL oder 125 g nicht mit H- und P-Sätzen gekennzeichnet werden.

Component	Hazard contents	GHS symbol	Hazard phrases	Precaution phrases
<i>Inhalt</i>	<i>Gefahrstoff</i>	<i>GHS-Symbol</i>	<i>H-Sätze</i>	<i>P-Sätze</i>
NucleoZOL	Phenol 30–60 % and guanidinium thiocyanate 30–60 % <i>Phenol 30–60 % und Guanidinthiocyanat 30–60 %</i> CAS 108-95-2, 593-84-0	 DANGER GEFAHR	301, 311, 314, 331, 341, 373, 412, EUH031	201, 202, 260, 273, 280, 301+310, 301+330+331, 302+352, 303+361+353, 304+340, 305+351+338, 308+313, 311, 361+364, 405, 501

Hazard phrases

H301	Toxic if swallowed. <i>Giftig bei Verschlucken.</i>
H311	Toxic in contact with skin. <i>Giftig bei Hautkontakt.</i>
H314	Causes severe skin burns and eye damage. <i>Verursacht schwere Verätzungen der Haut und schwere Augenschäden.</i>
H331	Toxic if inhaled. <i>Giftig bei Einatmen.</i>
H341	Suspected of causing genetic defects. <i>Kann vermutlich genetische Defekte verursachen.</i>
H373	May cause damage to organs through prolonged or repeated exposure. <i>Kann die Organe schädigen bei längerer oder wiederholter Exposition.</i>
H 412	Harmful to aquatic life with long lasting effects. <i>Schädlich für Wasserorganismen, mit langfristiger Wirkung.</i>
EUH031	Contact with acids liberates toxic gas. <i>Entwickelt bei Berührung mit Säure giftige Gase.</i>

Precaution phrases

- P201 Obtain special instructions before use.
Vor Gebrauch besondere Anweisungen einholen.
- P202 Do not handle until all safety precautions have been read and understood.
Vor Gebrauch alle Sicherheitshinweise lesen und verstehen.
- P260 Do not breathe vapors.
Dampf nicht einatmen.
- P273 Avoid release to the environment.
Freisetzung in die Umwelt vermeiden
- P280 Wear protective gloves / eye protection.
Schutzhandschuhe / Augenschutz tragen.
- P301+310 IF SWALLOWED: Immediately call a POISON CENTER / doctor / ...
BEI VERSCHLUCKEN: Sofort GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.
- P301+330+331 IF SWALLOWED: Rinse mouth. Do NOT induce vomiting.
BEI VERSCHLUCKEN: Mund ausspülen. KEIN Erbrechen herbeiführen.
- P302+352 IF ON SKIN: Wash with plenty of water / ...
BEI BERÜHRUNG MIT DER HAUT: Mit viel Wasser/... waschen.
- P303+361+353 IF ON SKIN (or hair): Take off immediately all contaminated clothing. Rinse skin with water / shower.
BEI BERÜHRUNG MIT DER HAUT (oder dem Haar): Alle kontaminierten Kleidungsstücke sofort ausziehen. Haut mit Wasser abwaschen / duschen.
- P304+340 IF INHALED: Remove person to fresh air and keep comfortable for breathing.
BEI EINATMEN: Die Person an die frische Luft bringen und für ungehinderte Atmung sorgen.
- P305+351+338 IF IN EYES: Rinse cautiously with water for several minutes.
BEI BERÜHRUNG MIT DEN AUGEN: Einige Minuten lang behutsam mit Wasser spülen.
- P308+313 IF exposed or concerned: Get medical advice / attention.
BEI Exposition oder falls betroffen: Ärztlichen Rat einholen / ärztliche Hilfe hinzuziehen.
- P311 Call a POISONCENTER / doctor / ...
GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.
- P361+364 Take off immediately all contaminated clothing and wash it before reuse.
Alle kontaminierten Kleidungsstücke sofort ausziehen und vor erneutem Tragen waschen.
- P405 Store locked up.
Unter Verschluss aufbewahren.
- P501 Dispose of contents / container to ...
Inhalt/Behälter ... zuführen.

For further information please see Material Safety Data Sheets (www.mn-net.com).
Weiterführende Informationen finden Sie in den Sicherheitsdatenblättern (www.mn-net.com).

5 NucleoZOL protocols

5.1 Isolation of small and large RNA in two separate fractions

Please note that RNA is separated in small RNA (10–200 nt) and large RNA (> 200 nt) in two fractions following this protocol.

1 Homogenize and lyse sample

Tissue

Homogenize tissue samples with a rotor-stator homogenizer or another mechanical disruption device using up to **100 mg of tissue per 1 mL NucleoZOL**. For tissues with high DNA content (e.g., spleen), it is recommended to use 50 mg of tissue/mL reagent.

For simplicity, this protocol describes RNA isolation using 1 mL of NucleoZOL. For processing of samples in 1.5 mL or 2 mL microcentrifuge tubes, use an 880 μ L aliquot of the homogenate (80 mg tissue + 800 μ L NucleoZOL). Residual homogenate can be stored at -20 °C or -70 °C for at least one year for later use.

Cells

Cells grown in monolayer: Remove cell culture medium and lyse cells by addition of at least **1 mL NucleoZOL** to the culture dish (diameter 3.5 cm, 10 cm²). Ensure complete lysis by repeated pipetting. Calculate the amount of reagent based on culture dish area, not on cell number.

An insufficient volume of the reagent will lead to DNA contamination of the isolated RNA.

Cells grown in suspension Sediment cells and lyse directly by the addition of NucleoZOL. Add at least **1 mL NucleoZOL per 10⁷ cells** and lyse cells by pipetting up and down several times.

Do not wash the cells before addition of NucleoZOL. Washing of cells might contribute to RNA degradation.

Liquid samples

Add **1 mL NucleoZOL per 400 μ L liquid sample** for homogenization and lysis. For processing sample volumes smaller than 400 μ L, add 1 mL of NucleoZOL and add RNase-free water to a final volume of 1.4 mL.

Fatty samples

Homogenize lipid-rich samples as described above. Centrifuge the samples for **5 min at 12,000 x g**. After centrifugation, a fat layer appears on top of the sample. Pierce the upper layer with a syringe/pipette tip and transfer the supernatant into a new tube.

2 Precipitate contaminants

Add **400 µL RNase-free water per 1 mL NucleoZOL** to the lysate. **Shake** the sample vigorously for **15 s**. **Incubate** at **room temperature (18–25 °C)** for **5–15 min**.

For samples containing 100 mg tissue/mL NucleoZOL, 15 min incubation at room temperature is recommended.

Centrifuge samples for **15 min** at **12,000 x g**.

Centrifugation can be performed at 4–28 °C.

A semi-solid pellet containing DNA, proteins, and polysaccharides forms at the bottom of the tube. The RNA is still solubilized in the supernatant.

3 Precipitate large RNA

Pipette **1 mL** of the supernatant to a new tube. Leave a layer of the supernatant on top of the precipitate. Add **400 µL 75 % ethanol** to **1 mL supernatant** for precipitation of the RNA.

Incubate samples at **room temperature** for **10 min**.

Centrifuge the samples for **8 min** at **12,000 x g**. A white pellet containing the RNA will be formed at the bottom of the tube. Transfer the supernatant containing the small RNA to a new tube and store it at 4 °C or at -20 °C.

The small RNA containing supernatant can be stored at -20 °C for one year.

4 Precipitate small RNA

Add **0.8 volumes of isopropanol** to the supernatant (~ 1.4 mL) obtained after precipitation of RNA (step 3) Incubate the samples for **30 min** at **4 °C**

Centrifuge the sample for **15 min** at **12,000 x g**.

Centrifugation can be performed at 4–28 °C.

Precipitated RNA will form a white pellet at the bottom of the tube.

5 Wash RNA**Large RNA**

Add **400–600 µL 75 % ethanol** when working with 1.5 mL tubes. For larger tubes, add **500 µL 75 % ethanol per 1 mL supernatant** used for precipitation.

Centrifuge for **1–3 min** at **4,000–8,000 x g**. Remove ethanol using a micropipette. Repeat washing step. **Do not dry the pellet.**

Small RNA

Add **400–600 µL 70 % isopropanol** when working with 1.5 mL tubes. For larger tubes, add **500 µL 70 % isopropanol per 1 mL supernatant** used for precipitation.

Centrifuge for **1–3 min** at **4,000–8,000 x g**. Remove ethanol using a micropipette. Repeat washing step. **Do not dry the pellet.**

Drying the RNA pellet may lead to a decrease in solubility.

6 Reconstitute RNA

Dissolve the RNA pellet in RNase-free water to obtain an RNA concentration of 1–2 µg/µL for the large RNA fraction and about 0.1 µg/µL for the small RNA fraction. Vortex the sample for **2–5 min** at **room temperature** for efficient solubilization.

For accurate determination of RNA concentration by OD measurement with a cuvette, dilute RNA in RNase-free water with a slightly alkaline pH, 1 mM NaOH or buffer with a pH > 8 (e.g., Elution Buffer AE, see ordering information 6.3). Distilled water typically has a pH < 7.

Note: The large RNA fraction contains RNA > 200 nt and contains 80–85 % of cellular RNA.

5.2 Isolation of total RNA

Total RNA (including small RNA, e.g., miRNA) is isolated with the following protocol.

1 Homogenization

Tissue

Homogenize tissue samples with a rotor-stator homogenizer or another mechanical disruption device using up to **100 mg of tissue per 1 mL NucleoZOL**. For tissues with high DNA content (e.g., spleen) it is recommended to use 50 mg of tissue/mL reagent.

For simplicity, this protocol describes RNA isolation using 1 mL of NucleoZOL. For processing the sample in 1.5 mL or 2 mL microcentrifuge tubes, use an 880 μ L aliquot of the homogenate (80 mg tissue + 800 μ L NucleoZOL). Residual homogenate can be stored at -20 °C or -70 °C for at least one year for later use.

Cells

Cells grown in monolayer: Remove cell culture medium and lyse cells by addition of at least **1 mL of NucleoZOL** to the culture disk (diameter 3.5 cm, 10 cm²). Ensure complete lysis by repeated pipetting. Calculate the amount of reagent based on culture dish area, not on cell number.

An insufficient volume of the reagent will lead to DNA contamination of the isolated RNA.

Cells grown in suspension: Sediment cells and lyse directly by the addition of NucleoZOL. Add at least **1 mL NucleoZOL per 10⁷ cells** and lyse cells by pipetting up and down several times.

Do not wash the cells before addition of NucleoZOL. Washing of cells might contribute to RNA degradation.

Liquid samples

Add **1 mL NucleoZOL per 400 μ L liquid sample** for homogenization and lysis. For processing sample volumes smaller than 400 μ L, add 1 mL of NucleoZOL and add water to a final volume of 1.4 mL.

Fatty samples

Homogenize lipid-rich samples as described above. Centrifuge the samples for **5 min at 12,000 x g**. After centrifugation, a fat layer appears on top of the sample. Pierce the upper layer with a syringe/pipette tip and transfer the supernatant into a new tube.

2 Precipitate contaminants

Add **400 µL RNase-free water per 1 mL NucleoZOL** to the lysate. **Shake** the sample vigorously for **15 s**. Incubate at **room temperature** for **5–15 min**.

For samples containing 100 mg tissue/mL NucleoZOL, 15 min incubation at room temperature is recommended.

Centrifuge samples for **15 min** at **12,000 x g**.

Centrifugation can be performed at 4–28 °C.

A semi-solid pellet containing DNA, proteins and polysaccharides forms at the bottom of the tube. The RNA is still solubilized in the supernatant.

Transfer **1 mL supernatant** (75% of total supernatant volume) to a fresh tube. Leave a layer of the supernatant above the DNA/protein pellet.

The pellet containing DNA, protein, and polysaccharides comprises approximately 10% in volume of the total homogenate-water mix (e.g., about 8% pellet for 80 mg tissue lysed in 1 mL reagent).

3 Phase separation (optional)

The basic protocol for total RNA isolation can be complemented by an optional phase separation. This is useful for samples with high DNA content and/or extracellular material.

Add **5 µL (0.5% of supernatant volume) 4-bromoanisole** to **1 mL transferred supernatant**. **Mix** well for 15 s and incubate at **room temperature** for **35 min**.

Do not substitute 4-bromoanisole with bromchloropropane or chloroform!

Centrifuge for **10 min** at **12,000 x g** (4–25 °C).

Residual DNA, proteins, and polysaccharides accumulate in the organic phase at the bottom of the tube. RNA is still solubilized in the supernatant.

4 Precipitate total RNA

Pipette RNA containing supernatant from step 2 or 3 into a fresh tube.

Add **1 mL of isopropanol** per **1 mL supernatant** in order to precipitate RNA. Incubate samples at **room temperature** for **10 min**.

Centrifuge samples for **10 min** at **12,000 x g**.

Typically, RNA is obtained as a white pellet at the bottom of the tube. For spleen samples, RNA forms a gel-like membrane on the bottom of the tube. Upon washing with ethanol, the membrane becomes more visible.

5 Wash RNA

Use **400–600 µL 75 % ethanol** when precipitating in 1.5 mL tubes. For larger tubes, add **500 µL 75 % ethanol per 1 mL supernatant**

Centrifuge the pellets **for 1–3 min at 4,000–8,000 x g**. Remove ethanol from the pellet by pipetting. Repeat the ethanol washing step. **Do not dry the RNA pellet between wash steps.**

Drying the RNA pellet may lead to a decrease in solubility.

6 Reconstitute RNA

Dissolve the RNA pellet in RNase-free water to obtain an RNA concentration of 1–2 µg/µL. Vortex the sample **2–5 min at room temperature** for efficient solubilization.

For accurate determination of RNA concentration by OD measurement with a cuvette, dilute RNA in RNase-free water with a slightly alkaline pH, 1 mM NaOH or buffer with a pH > 8 (e.g., Elution Buffer AE, see ordering information, section 6.3). Distilled water typically has a pH < 7.

5.3 Isolation of total RNA in combination with NucleoSpin® RNA kits

1 Homogenization

The homogenization is similar to the standard protocols. For detailed information refer to section 5.1 or 5.2.

Tissue

Use up to **50 mg** tissue with **500 µL NucleoZOL**. Larger samples may exceed the RNA binding capacity of the NucleoSpin® RNA Binding Column.

Cells

Use up to **10⁷ cells** with **1 mL NucleoZOL** per preparation.

Liquid samples

Use up to **100 µL liquid sample** with **250 µL NucleoZOL**.

2 Precipitate contaminants

Add **400 µL RNase-free water** to the lysate per **1 mL NucleoZOL** used for homogenization/lysis.

Shake the sample vigorously for **15 s** and incubate at **room temperature** for **5–15 min**.

For samples containing 100 mg tissue/mL NucleoZOL, 15 min incubation at room temperature is recommended.

Centrifuge samples for **15 min** at **12,000 x g**.

Centrifugation can be performed at 4–28 °C.

A semi-solid pellet containing DNA, proteins, and polysaccharides forms at the bottom of the tube. The RNA is still solubilized in the supernatant.

Transfer **500 µL supernatant** into a fresh tube. Do not disturb the DNA/protein pellet.

The pellet containing DNA, protein, and polysaccharides comprises approximately 10% in volume of the total homogenate-water mix (e.g., about 8% pellet for 80 mg tissue lyses in 1 mL reagent).

3 Adjust RNA binding conditions

Add **1 vol 70% ethanol** to **1 vol NucleoZOL** sample homogenate and mix by vortexing.

4 Bind RNA

For each preparation, take one **NucleoSpin® RNA Column** (light blue ring) placed in a Collection Tube and load the lysate. Centrifuge for **30 s** at **8,000 x g**.

Discard Collection Tube with flow-through and place the column in a new Collection Tube.

Maximum loading capacity of NucleoSpin® RNA Columns is 750 µL. Repeat the procedure if larger volumes are processed.

5 Wash and dry silica membrane

1st wash

Add **700 µL Buffer RA3** to the NucleoSpin® RNA Column. Centrifuge for **30 s** at **8,000 x g**. Discard flow-through and reuse Collection Tube.

2nd wash

Add **350 µL Buffer RA3** to the NucleoSpin® RNA Column. Centrifuge for **2 min** at **8,000 x g**.

Place the NucleoSpin® RNA Column into an RNase-free Collection Tube (1.5 mL, supplied). Open the lid of the column and let the membrane **dry** for **3 min**.

If the liquid level in the Collection Tube has reached the NucleoSpin® RNA Column after centrifugation, discard flow-through and centrifuge again.

The procedure ensures complete removal of ethanol from the column.

6 Elute RNA

Add **60 µL RNase-free H₂O** onto the center of the membrane and centrifuge for **1 min** at **8,000 x g**.

6 Appendix

6.1 Digestion of residual DNA solution

NucleoZOL efficiently removes DNA when processing samples according to the standard protocol, resulting in minimal residual DNA in the purified RNA. Residual DNA will not be detectable in most downstream applications. If large samples or samples with high levels of DNA are processed, it may be difficult to remove all traces of DNA. The amount of residual DNA depends on the sample type, amount, DNA content and the detection sensitivity of the method used to analyze residual DNA. A typical example is an RT-PCR reaction in which the primer molecules do not differentiate between cDNA (derived from RNA) and contaminating genomic DNA. The effect is prominent if

- high copy number targets are analyzed (e.g., multi gene family, mitochondrial, plastid or plasmid targets (from transfections)
- the target gene is of a very low expression level
- the amplicon is relatively small (< 200 bp).

DNA digestion in solution can efficiently destroy contaminating DNA. However, stringent RNase control and subsequent repurification of the RNA (in order to remove buffer, salts, DNase and digested DNA) are usually required. High quality, RNase-free, recombinant rDNase (REF 740963, see ordering information 6.3) facilitates such a digestion in solution in order to remove traces of contaminating DNA.

A) Digest DNA (Reaction setup)

Add **6 µL Reaction Buffer for rDNase** and **0.6 µL rDNase** to **60 µL eluted RNA**. (Alternatively, premix 100 µL Reaction Buffer for rDNase and 10 µL rDNase and add 1/10 volume to one volume of RNA eluate). Gently swirl the tube in order to mix the solution. Spin down gently (approx. 1 s at 1,000 x g) to collect every droplet of the solution at the bottom of the tube.

B) Incubate sample

Incubate for **10 min** at **37 °C**.

C) Repurify RNA

Repurify RNA with a suitable RNA cleanup procedure: NucleoSpin® RNA Clean-up, NucleoSpin® RNA Clean-up XS kits (see ordering information 6.3), or by ethanol precipitation.

Ethanol precipitation, exemplary:

Add **0.1 volume** of **3 M sodium acetate, pH 5.2** and **2.5 volumes** of **96–100 % ethanol** to **one volume of sample**. Mix thoroughly. Incubate **several minutes** to **several hours** at **-20 °C** or **4 °C**.

Note: Choose long incubation times if the sample contains low RNA concentration. Short incubation times are sufficient if the sample contains high RNA concentration.

Centrifuge for **10 min** at **maximum speed**.

Wash RNA pellet with **70 % ethanol**.

Dry RNA pellet and resuspend RNA in RNase-free water.

6.2 Troubleshooting

Problem	Possible cause and suggestions
Insufficient yield	<i>Homogenization or sample lysis is incomplete</i>
	<ul style="list-style-type: none"> • Improve homogenization by testing more stringent conditions.
	<i>Solubilization of the RNA pellet is incomplete.</i>
	<ul style="list-style-type: none"> • Increase volume of RNase-free water for dissolving of RNA and prolong mixing for solubilization.
Ratio $A_{260/280}$ < 1.6	<i>Volume of NucleoZOL used for homogenization was too low</i>
	<ul style="list-style-type: none"> • Increase volume of NucleoZOL.
	<i>Low pH during spectrophotometric quantification</i>
	<ul style="list-style-type: none"> • Use Buffer AE (5 mM Tris pH 8.5) for sample dilution for spectrophotometric quantification.
	<i>RNA pellet was only partly solubilized</i>
	<ul style="list-style-type: none"> • Increase volume of RNase-free water for dissolving of RNA and prolong mixing for solubilization.
	<i>Contamination of polysaccharide or proteoglycan</i>
	<ul style="list-style-type: none"> • Perform phase separation as described in section 5.2.3.
Degraded RNA	<i>Inadequate tissue sampling</i>
	<ul style="list-style-type: none"> • Make sure to use fresh tissue or flash-freeze tissue immediately upon harvest.
	<i>Inappropriate storage conditions</i>
	<ul style="list-style-type: none"> • Store samples at -70 °C.
	<i>Cell dissolution during trypsinization</i>
	<ul style="list-style-type: none"> • Make sure cells stay intact during trypsinization.
	<i>RNase contaminated solutions or tubes</i>
	<ul style="list-style-type: none"> • Make sure to work in an RNase-free environment.

Problem	Possible cause and suggestions
Contamination with DNA	<p><i>Volume of NucleoZOL used for homogenization was too low</i></p> <ul style="list-style-type: none"> • Increase volume of NucleoZOL.
	<p><i>Sample material contains strong buffers, organic solvents, alkaline solution, or salt</i></p> <ul style="list-style-type: none"> • The precipitation of DNA (step 2) can be improved by the following modification: Increase incubation time to 15 min after addition of water (step 2). Centrifuge at 16,000 x g. • Use support protocol 6.1 for subsequent rDNase digestion in solution
Contamination with proteoglycan, fat, or polysaccharide	<p><i>Inefficient precipitation of contaminants</i></p> <ul style="list-style-type: none"> • Centrifuge the initial crude homogenate (step 1) for phase separation in an additional step for 10 min at 12,000 x g.

6.3 Ordering information

Product	REF	Pack of
NucleoZOL	740404.200	200 mL
NucleoSpin® RNA	740955.10/50/250	10/50/250 preps
NucleoSpin® RNA Plus	740984.10/50/250	10/50/250 preps
NucleoSpin® RNA XS	740902.10/50/250	10/50/250 preps
NucleoSpin® RNA Clean-up XS	740903.10/50/250	10/50/250 preps
NucleoSpin® RNA Clean-up	740948.10/50/250	10/50/250 preps
RNase-free Water	740378.1000	1000 mL
Elution Buffer AE	740917.1	1000 mL
rDNase Set	740963	one set

6.4 Product use restriction / warranty

NucleoZOL kit components are intended, developed, designed, and sold FOR RESEARCH PURPOSES ONLY, except, however, any other function of the product being expressly described in original MACHEREY-NAGEL product leaflets.

MACHEREY-NAGEL products are intended for GENERAL LABORATORY USE ONLY! MACHEREY-NAGEL products are suited for QUALIFIED PERSONNEL ONLY! MACHEREY-NAGEL products shall in any event only be used wearing adequate PROTECTIVE CLOTHING. For detailed information please refer to the respective Material Safety Data Sheet of the product! MACHEREY-NAGEL products shall exclusively be used in an ADEQUATE TEST ENVIRONMENT. MACHEREY-NAGEL does not assume any responsibility for damages due to improper application of our products in other fields of application. Application on the human body is STRICTLY FORBIDDEN. The respective user is liable for any and all damages resulting from such application.

DNA/RNA/PROTEIN purification products of MACHEREY-NAGEL are suitable for IN-VITRO-USES ONLY!

ONLY MACHEREY-NAGEL products specially labeled as IVD are also suitable for IN-VITRO-diagnostic use. Please pay attention to the package of the product. IN-VITRO-diagnostic products are expressly marked as IVD on the packaging.

IF THERE IS NO IVD SIGN, THE PRODUCT SHALL NOT BE SUITABLE FOR IN-VITRO-DIAGNOSTIC USE!

ALL OTHER PRODUCTS NOT LABELED AS IVD ARE NOT SUITED FOR ANY CLINICAL USE (INCLUDING, BUT NOT LIMITED TO DIAGNOSTIC, THERAPEUTIC AND/OR PROGNOSTIC USE).

No claim or representations is intended for its use to identify any specific organism or for clinical use (included, but not limited to diagnostic, prognostic, therapeutic, or blood banking). It is rather in the responsibility of the user or - in any case of resale of the products - in the responsibility of the reseller to inspect and assure the use of the DNA/RNA/protein purification products of MACHEREY-NAGEL for a well-defined and specific application.

MACHEREY-NAGEL shall only be responsible for the product specifications and the performance range of MN products according to the specifications of in-house quality control, product documentation and marketing material.

This MACHEREY-NAGEL product is shipped with documentation stating specifications and other technical information. MACHEREY-NAGEL warrants to meet the stated specifications. MACHEREY-NAGEL's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Supplementary reference is made to the general business terms and conditions of MACHEREY-NAGEL, which are printed on the price list. Please contact us if you wish to get an extra copy.

There is no warranty for and MACHEREY-NAGEL is not liable for damages or defects arising in shipping and handling (transport insurance for customers excluded), or out of accident or improper or abnormal use of this product; defects in products or

components not manufactured by MACHEREY-NAGEL, or damages resulting from such non-MACHEREY-NAGEL components or products.

MACHEREY-NAGEL makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, REPRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO MACHEREY-NAGEL PRODUCTS.

In no event shall MACHEREY-NAGEL be liable for claims for any other damages, whether direct, indirect, incidental, compensatory, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of MACHEREY-NAGEL products to perform in accordance with the stated specifications. This warranty is exclusive and MACHEREY-NAGEL makes no other warranty expressed or implied.

The warranty provided herein and the data, specifications and descriptions of this MACHEREY-NAGEL product appearing in MACHEREY-NAGEL published catalogues and product literature are MACHEREY-NAGEL's sole representations concerning the product and warranty. No other statements or representations, written or oral, by MACHEREY-NAGEL's employees, agent or representatives, except written statements signed by a duly authorized officer of MACHEREY-NAGEL are authorized; they should not be relied upon by the customer and are not a part of the contract of sale or of this warranty.

Product claims are subject to change. Therefore please contact our Technical Service Team for the most up-to-date information on MACHEREY-NAGEL products. You may also contact your local distributor for general scientific information. Applications mentioned in MACHEREY-NAGEL literature are provided for informational purposes only. MACHEREY-NAGEL does not warrant that all applications have been tested in MACHEREY-NAGEL laboratories using MACHEREY-NAGEL products. MACHEREY-NAGEL does not warrant the correctness of any of those applications.

Last updated: 07/2010, Rev. 03

Please contact:
MACHEREY-NAGEL GmbH & Co. KG
Tel.: +49 24 21 969-270
tech-bio@mn-net.com

Trademarks/patents/disclaimer:

NucleoSpin® is a registered trademark of MACHEREY-NAGEL GmbH & Co KG
RNAlater® is a trademark of Ambion Inc.

NucleoZOL reagent: US Patents 7,794,932 and 8,367,817 B2 and international patents

All used names and denotations can be brands, trademarks, or registered labels of their respective owner – also if they are not special denotation. To mention products and brands is only a kind of information (i.e., it does not offend against trademarks and brands and can not be seen as a kind of recommendation or assessment). Regarding these products or services we can not grant any guarantees regarding selection, efficiency, or operation.