



# RNA from blood

## User manual

NucleoSpin<sup>®</sup> RNA Blood  
NucleoSpin<sup>®</sup> RNA Blood Midi

This product distributed by  
**Takara Bio USA, Inc.**

**Ordering information**  
1.800.662.2566  
ordersUS@takarabio.com

**Technical support**  
1.800.662.2566  
techUS@takarabio.com



Clontech Takara cellartis













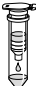





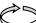
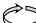



[takarabio.com](http://takarabio.com)

For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Takara Bio Inc. ©2018

September 2017 / Rev. 05

# RNA from blood

## Protocol-at-a-glance (Rev. 05)

	Mini NucleoSpin® RNA Blood 200 µL blood	Mini NucleoSpin® RNA Blood 400 µL blood	Midi NucleoSpin® RNA Blood Midi 1.3 mL blood
<b>1 Lyse blood</b>	 200 µL blood 200 µL DL Mix 5 µL Pro. K RT, 3–15 min (shaking)	 400 µL blood 400 µL DL Mix 10 µL Pro. K RT, 3–15 min (shaking)	 1.3 mL blood 1.3 mL DL Mix 33 µL Pro. K RT, 3–15 min (shaking)
<b>2 Adjust RNA binding conditions</b>	200 µL 70% ethanol Mix	400 µL 70% ethanol Mix	1.3 mL 70% ethanol Mix
<b>3 Bind RNA</b>	 Load sample 11,000 x g, 30 s	 Load sample stepwise 11,000 x g, 30 s	 Load sample 4,500 x g, 3 min
<b>4 Desalt silica membrane</b>	 350 µL MDB 11,000 x g, 30 s	 350 µL MDB 11,000 x g, 30 s	 1.2 mL MDB 4,500 x g, 3 min
<b>5 Digest DNA</b>	 95 µL rDNase RT, 15 min	 95 µL rDNase RT, 15 min	 240 µL rDNase RT, 15 min
<b>6 Wash silica membrane</b>	 200 µL RB2 600 µL RB3 250 µL RB3	 200 µL RB2 600 µL RB3 250 µL RB3	 1 mL RB2 3 mL RB3
<b>1<sup>st</sup> and 2<sup>nd</sup> wash</b>	 11,000 x g, 30 s	 11,000 x g, 30 s	 4,500 x g, 3 min
<b>3<sup>rd</sup> wash</b>	 11,000 x g, 2 min	 11,000 x g, 2 min	
<b>7 Elute RNA</b>	 60 µL RNase-free H <sub>2</sub> O 11,000 x g, 30 s	 60 µL RNase-free H <sub>2</sub> O 11,000 x g, 30 s	 200 µL RNase-free H <sub>2</sub> O 4,500 x g, 3 min

---

## Table of contents

1	Components	4
1.1	Kit contents	4
1.2	Reagents, consumables, and equipment to be supplied by user	6
1.3	About this user manual	6
2	Product description	7
2.1	The basic principle	7
2.2	Kit specifications	8
2.3	Handling, preparation, and storage of starting materials	9
2.4	Elution procedures	9
3	Storage conditions and preparation of working solutions	10
4	Safety instructions	11
5	NucleoSpin® RNA Blood protocols	13
5.1	RNA isolation from 200 µL blood	13
5.2	RNA isolation from 400 µL blood	16
6	NucleoSpin® RNA Blood Midi protocol – RNA isolation from 1.3 mL blood	19
7	Appendix	22
7.1	rDNase digestion in solution	22
7.2	Troubleshooting	24
7.3	Ordering information	27
7.4	Product use restriction/warranty	27

# 1 Components

## 1.1 Kit contents

<b>NucleoSpin® RNA Blood*</b>		
<b>REF</b>	<b>10 preps 740200.10</b>	<b>50 preps 740200.50</b>
Lysis Buffer DL	25 mL	25 mL
Wash Buffer RB2	13 mL	13 mL
Wash Buffer RB3 (Concentrate)**	6 mL	12 mL
Membrane Desalting Buffer MDB	10 mL	25 mL
Reaction Buffer for rDNase	7 mL	7 mL
rDNase, RNase-free (lyophilized)**	1 vial (size C)	2 vials (size D)
Liquid Proteinase K	120 µL	600 µL
RNase-free H <sub>2</sub> O	13 mL	13 mL
NucleoSpin® RNA Blood Columns (light blue rings - plus Collection Tubes)	10	50
Collection Tubes (2 mL, with lid) for lysis	10	50
Collection Tubes (1.5 mL) for elution	10	50
Collection Tubes (2 mL)	30	150
User manual	1	1

---

\* Patent pending

\*\* For preparation of working solutions and storage conditions see section 3.

## Kit contents continued

<b>NucleoSpin® RNA Blood Midi*</b>	
<b>REF</b>	<b>20 preps 740210.20</b>
Lysis Buffer DL	50 mL
Wash Buffer RB2	2 x 13 mL
Wash Buffer RB3 (Concentrate)**	25 mL
Membrane Desalting Buffer MDB	50 mL
Reaction Buffer for rDNase	7 mL
rDNase, RNase-free (lyophilized)**	2 vials (size D)
Liquid Proteinase K	800 µL
RNase-free H <sub>2</sub> O	13 mL
NucleoSpin® RNA Blood Midi Columns (plus Collection Tubes)	20
Collection Tubes (15 mL) for lysis, elution, and washing steps	60
User manual	1

\* Patent pending

\*\* For preparation of working solutions and storage conditions see section 3.

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

### Reagents

- 96–100 % ethanol (to prepare Wash Buffer RB3)
- 70 % ethanol (to adjust RNA binding conditions)

### Consumables

- Sterile RNase-free tips

### Equipment

- Manual pipettors
- Vortex mixer
- Centrifuge for microcentrifuge tubes (NucleoSpin® RNA Blood)
- Centrifuge for 15 mL tubes with a swing-out rotor capable of reaching 4,500 x *g* (NucleoSpin® RNA Blood Midi)
- Personal protection equipment (e.g., lab coat, gloves, goggles)

## 1.3 About this user manual

It is strongly recommended to read the detailed protocol sections of this user manual if using the **NucleoSpin® RNA Blood** or **NucleoSpin® RNA Blood Midi** kits for the first time. However, experienced users may refer to the Protocol-at-a-glance. The Protocol-at-a-glance is designed to be used only as a supplemental tool for quick referencing while performing the purification procedure.

All technical literature is available on the internet at [www.mn-net.com](http://www.mn-net.com).

## 2 Product description

### 2.1 The basic principle

The **NucleoSpin® RNA Blood** kits offer a direct total blood lysis from 200–400 µL (NucleoSpin® RNA Blood) or 400–1300 µL (NucleoSpin® RNA Blood Midi) whole blood collected in standard (e.g., EDTA) blood collection tubes. One of the most important aspects in RNA purification is to prevent RNA degradation during the isolation. With the **NucleoSpin® RNA Blood** method, leukocytes (the main source of RNA in whole blood) and other blood cells, are lysed by incubating the whole blood in a solution containing large amounts of chaotropic ions. This lysis buffer immediately inactivates RNases (which are present in virtually all biological materials) and creates appropriate binding conditions that favor adsorption of RNA to the silica membrane. A complex selective erythrocyte lysis and preparation of a leukocyte pellet is not necessary. Contaminating DNA, which is also bound to the silica membrane, is removed by a recombinant DNase solution (supplied) which is directly applied onto the silica membrane during the preparation. Simple washing steps with two different buffers remove salts, metabolites, and macromolecular cellular components. Pure RNA is finally eluted under low ionic strength conditions with RNase-free H<sub>2</sub>O (supplied).

The RNA preparation using **NucleoSpin® RNA Blood** kits is performed at room temperature. A refrigerated centrifuge is not necessary. The eluate, however, should be handled with care because RNA is very sensitive to trace contaminations of RNases, often found on general lab ware, fingerprints, and dust. To ensure RNA stability, keep RNA frozen at -20 °C for short-term or -70 °C for long-term storage.

#### **Simultaneous isolation of RNA and DNA (NucleoSpin® RNA/DNA Buffer Set)**

The NucleoSpin® RNA/DNA Buffer Set (see ordering information) is a support set for RNA and DNA isolation in conjunction with NucleoSpin® RNA, NucleoSpin® RNA XS, NucleoSpin® miRNA, NucleoSpin® RNA Plant, NucleoSpin® RNA/Protein, and NucleoSpin® RNA Blood.

This patented technology enables successive elution of DNA and RNA from one NucleoSpin® Column with low salt buffer and water respectively. DNA and RNA are immediately ready for downstream applications.

## 2.2 Kit specifications

- **NucleoSpin® RNA Blood** kits are recommended for the isolation of RNA from whole blood (e.g., stabilized with EDTA, citrate, or heparin).
- The **NucleoSpin® RNA Blood** kits allow the purification of RNA with an  $A_{260}/A_{280}$  ratio typically exceeding 1.9 (measured in TE buffer, pH 7.5).
- The isolated RNA is ready to use for typical downstream applications (e.g., reverse transcriptase-PCR (RT-PCR)).
- RNA isolated with the NucleoSpin® RNA Blood kits is typically of high integrity. However, RNA integrity strongly depends on the sample quality.
- The amount of DNA contamination is significantly reduced during on-column digestion with rDNase. However, in very sensitive applications, it may be possible to detect traces of DNA. The probability of DNA detection with PCR increases with:
  1. the number of DNA copies per preparation: single copy target < plasmidial / mitochondrial target < plasmid transfected into cells.
  2. decreasing PCR amplicon size.

**Table 1: Kit specifications at a glance**

Parameter	NucleoSpin® RNA Blood	NucleoSpin® RNA Blood Midi
Sample material	200–400 µL fresh or frozen whole blood (e.g., stabilized with EDTA, citrate, or heparin)	400-1300 µL fresh or frozen whole blood (e.g., stabilized with EDTA, citrate, or heparin)
Format	Mini spin column	Midi spin column
Fragment size	> 200 nt	> 200 nt
Typical yield	~ 7 µg (3–20 µg) per 1 mL blood from healthy subjects	~ 7 µg (3–20 µg) per 1 mL blood from healthy subjects
$A_{260}/A_{280}$	1.9–2.1	1.9–2.1
Elution volume	40–120 µL	200–400 µL
Binding capacity	200 µg	700 µg
Preparation time	45 min/6 preps (excl. lysis)	85 min/6 preps (excl. lysis)

The **NucleoSpin® RNA Blood** kit contains one protocol that allows the use of 200 µL of whole blood by a total direct blood lysis and a second protocol for processing 400 µL of whole blood with a second loading step.

The NucleoSpin® RNA Blood Midi kit contains a protocol that allows 1.3 mL of whole blood by a total direct blood lysis.

If other volumes than 200 µL, 400 µL, or 1300 µL blood are used, adjust the volumes of Buffer DL and 70 % ethanol in step 1 and 2 of the corresponding protocol by maintaining the following ratio:

**1:1:1** (sample / Buffer DL / 70 % ethanol)

Example: 300 µL blood + 300 µL Buffer DL + 300 µL 70 % ethanol



The volume of Proteinase K can be calculated as follows:

Blood volume  $\mu\text{L}$  / 40 = volume Proteinase K  $\mu\text{L}$

**Example:** 300  $\mu\text{L}$  blood / 40 = 7.5  $\mu\text{L}$  Liquid Proteinase K

The isolated RNA can be used as a template in RT-PCR-reactions. Generally, 1–40% of the eluate from RNA prepared with 200–400  $\mu\text{L}$  blood is suitable as a template for RT-PCR. If possible, intron-spanning primers should be used for RT-PCR.

## 2.3 Handling, preparation, and storage of starting materials

**NucleoSpin® RNA Blood** kits are designed for isolation of total RNA from fresh, human whole blood. Whole blood should be collected in the presence of an anticoagulant, preferably EDTA, citrate, or heparin.

It is highly recommended to process blood samples within a few hours after collecting them (when EDTA, citrate, or heparin collection tubes are used). Samples should be stored at 4 °C for no longer than 24 hours. The mRNAs contained in blood cells have different stabilities. As a result, in order to ensure that the isolated RNA contains a representative distribution of mRNAs, blood samples should not be stored for long periods before isolating RNA.

If frozen blood samples have to be processed, aliquots of 200  $\mu\text{L}$ , 400  $\mu\text{L}$ , or 1300  $\mu\text{L}$ , preferably, of frozen blood aliquots should be quickly thawed in the presence of 1 volume Lysis Buffer DL while shaking.

If long-term storage of stabilized whole blood is necessary, it is recommended storing the lysates at -20 °C. For this, add the indicated volume of Lysis Buffer DL to the blood sample without adding Liquid Proteinase K. Store the lysates at -20 °C. After thawing, add Liquid Proteinase K and follow the protocol at step 1.

Wear gloves at all times during the preparation. Change gloves frequently.

## 2.4 Elution procedures

It is possible to adjust the elution method and volume of RNase-free water used for the subsequent application of interest (refer to Table 1 regarding suitable ranges of elution volumes). In addition to the standard method described in the individual protocols (recovery rate about 70–90%), there are several modifications possible:

- **High yield:** Perform two elution steps with the volume indicated in the individual protocol. About 90–100% of bound nucleic acids will be eluted.
- **High yield and high concentration:** Elute with the standard elution volume and apply the eluate once more onto the column for re-elution.

Eluted RNA should be immediately placed and kept on ice for optimal stability and to prohibit omnipresent RNases (general lab ware, fingerprints, dust) from degrading the RNA. For short term storage, freeze at -20 °C, for long term storage, freeze at -70 °C.

### 3 Storage conditions and preparation of working solutions

**Attention:** Buffers DL, RB2, and MDB contain chaotropic salts. Wear gloves and goggles!

**CAUTION:** Buffers DL, RB2, and MDB contain guanidinium salts which can form highly reactive compounds when combined with bleach (sodium hypochlorite). DO NOT add bleach or acidic solutions directly to the sample-preparation waste.

- Store lyophilized **rDNase (RNase-free)** at 4 °C on arrival (stable up to 1 year).
- Lysis Buffer DL is light sensitive during long time storage. Therefore, Lysis Buffer DL is provided in a black bottle. Short light exposure (several hours) does not affect the buffer.
- After first use, it is recommended to store Liquid Proteinase K at 4 °C or -20 °C.
- All other kit components should be stored at room temperature (18–25 °C) and are stable up to one year. Storage at lower temperatures may cause precipitation of salts.
- Check that 70 % ethanol is available as additional solution to adjust RNA binding conditions.
- Check that 96–100 % ethanol is available as additional solution to prepare Wash Buffer RB3.

Before starting any **NucleoSpin® RNA Blood** protocol, prepare the following:

- **rDNase (RNase-free):** Add indicated volume of Reaction Buffer for rDNase (see table below) to the rDNase vial and incubate for 1 min at room temperature. Gently swirl the vials to completely dissolve the rDNase. Be careful not to mix rDNase vigorously as rDNase is sensitive to mechanical agitation. Dispense into aliquots and store at -20 °C. The frozen working solution is stable for 6 months. Do not freeze/thaw the aliquots more than three times. (Be careful when opening the vial as some particles of the lyophilisate may be attached to the lid.)
- **Wash Buffer RB3:** Add the indicated volume of 96–100 % ethanol (see table below) to Buffer RB3 Concentrate. Mark the label of the bottle to indicate that ethanol was added. Store Wash Buffer RB3 at room temperature (18–25 °C) for up to one year.

REF	NucleoSpin® RNA Blood		NucleoSpin® RNA Blood Midi
	10 preps 740200.10	50 preps 740200.50	20 preps 740210.20
Wash Buffer RB3 Concentrate	6 mL Add 24 mL ethanol	12 mL Add 48 mL ethanol	25 mL Add 100 mL ethanol
rDNase, RNase-free (lyophilized)	1 vial (size C) Add 1 mL Reaction Buffer for rDNase	2 vials (size D) Add 2.5 mL Reaction Buffer for rDNase to each vial	2 vials (size D) Add 2.5 mL Reaction Buffer for rDNase to each vial






## 4 Safety instructions

The following components of the **NucleoSpin® RNA Blood** and **NucleoSpin® RNA Blood Midi** kits contain 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>
DL	Guanidinium thiocyanate 30–60 % <i>Guanidinthiocyanat 30–60 %</i>  CAS 593-84-0	 <i>WARNING</i> <i>ACHTUNG</i>	302, 412, EUH031	260, 273
Liquid Proteinase K	Proteinase K 1–3 % <i>Proteinase K 1–3 %</i>  CAS 39450-01-6	 <i>WARNING</i> <i>ACHTUNG</i>	317	261, 280
MDB	Ethanol 5–20 % <i>Ethanol 5–20 %</i>  CAS 64-17-5	 <i>WARNING</i> <i>ACHTUNG</i>	226	210
RB2	Guanidine hydrochloride 24–36 % and ethanol 20–35 % <i>Guanidinhydrochlorid 24–36 %</i> <i>und Ethanol 20–35 %</i>  CAS 50-01-1, 64-17-5	 <i>WARNING</i> <i>ACHTUNG</i>	226, 302	210
rDNase	rDNase 90–100 % <i>rDNase 90–100 %</i>  CAS 9003-98-9	 <i>DANGER</i> <i>GEFAHR</i>	317, 334	261, 280, 342+311

### Hazard phrases

H 226	Flammable liquid and vapour. <i>Flüssigkeit und Dampf entzündbar.</i>
H 302	Harmful if swallowed. <i>Gesundheitsschädlich bei Verschlucken.</i>
H 317	May cause an allergic skin reaction. <i>Kann allergische Hautreaktionen verursachen.</i>
H 334	May cause allergy or asthma symptoms or breathing difficulties if inhaled. <i>Kann bei Einatmen Allergie, asthmaartige Symptome oder Atembeschwerden verursachen.</i>

H 412 Harmful to aquatic life with long lasting effects.  
*Schädlich für Wasserorganismen, mit langfristiger Wirkung.*

EUH 031 Contact with acids liberates toxic gas.  
*Entwickelt bei Berührung mit Säure giftige Gase.*

### Precaution phrases

P 210 Keep away from heat/sparks/open flames/hot surfaces.  
*Von Hitze, heißen Oberflächen, Funken, offenen Flammen sowie anderen Zündquellenarten fernhalten.*

P 260 In contact with water releases flammable gases which may ignite spontaneously.  
*In Berührung mit Wasser entstehen entzündbare Gase, die sich spontan entzünden können.*

P 261 In contact with water releases flammable gases.  
*In Berührung mit Wasser entstehen entzündbare Gase.*

P 273 Avoid release to the environment.  
*Freisetzung in die Umwelt vermeiden.*

P 280 Wear protective gloves / protective clothing / eye protection / face protection.  
*Schutzhandschuhe / Schutzkleidung / Augenschutz / Gesichtsschutz tragen.*

P 342 If experiencing respiratory symptoms:  
*Bei Symptomen der Atemwege:*

P 342+311 If experiencing respiratory symptoms: Call a POISON CENTER / doctor / ...  
*Bei Symptomen der Atemwege: GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.*



The symbol shown on labels refers to further safety information in this section.  
*Das auf Etiketten dargestellte Symbol weist auf weitere Sicherheitsinformationen dieses Kapitels hin.*

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

## 5 NucleoSpin® RNA Blood protocols

### 5.1 RNA isolation from 200 µL blood

Before starting the preparation:

- Check if Wash Buffer RB3 and rDNase were prepared according to section 3.
- The complete procedure should be performed at room temperature (18–25 °C).
- The use of blood collected in common blood collection tubes with anticoagulant (typically EDTA) is recommended. For frozen blood samples, see section 2.3.
- Check if 70 % ethanol is available to adjust binding conditions.
- Refer to section 2.2 if less than 200 µL whole blood is used.

#### 1 Lyse blood

Provide **200 µL whole blood** in a Collection Tube (2 mL, with lid; provided).



**200 µL blood**

Add **200 µL Lysis Buffer DL** to the tube and close the lid. Mix. If necessary, shortly spin to clean the lid.

**+ 200 µL DL**

Add **5 µL Liquid Proteinase K** and close the lid.

**+ 5 µL Liquid Proteinase K**

Incubate **3–15 min** at **room temperature** (18–25 °C) vigorously shaking the tube on a shaker (e.g., Eppendorf Thermoshaker, 1,400 rpm) .

**RT, 3–15 min**

Centrifuge briefly to clean the lid (~ 1 s at ~ **2,000 x g**). Short spin only!



**~ 2,000 x g,  
~ 1 s**

#### 2 Adjust RNA binding conditions

Add **200 µL 70 % ethanol** to the tube and mix vigorously.



**+ 200 µL  
70 % ethanol**

*Note: It is important to thoroughly mix the ethanol with the lysate. Recommended: Place tubes in a rack with lid. Close the rack lid and strongly shake the assembly. Alternatively, pipette the solution up and down ~ 5 times.*

**Mix**

Centrifuge briefly to clean the lid (~ 1 s at ~ **2,000 x g**). Short spin only!



**~ 2,000 x g,  
~ 1 s**

**3 Bind RNA**

Adjust pipette to **610 µL** and transfer lysate into a **NucleoSpin® RNA Blood Column** placed in a Collection Tube.



**Load lysate**

*Note: Do not pipette more than 650 µL into the spin column, this will cause the column to overflow! Avoid formation of foam and aerosols! Avoid wetting the rim (edge) of the column.*

Centrifuge **30 s** at **11,000 x g**. Discard flow-through and Collection Tube. Place the column in a new Collection Tube (2 mL; provided).



**11,000 x g,  
30 s**

**4 Desalt silica membrane**

Add **350 µL MDB** (Membrane Desalting Buffer) onto the column and centrifuge **30 s** at **11,000 x g**.



**+ 350 µL MDB**

*Note: After centrifugation, the column can remain in the Collection Tube including the flow-through! The flow-through may be slightly brown. The flow-through can remain in the tube without disturbing DNA digestion.*



**11,000 x g,  
30 s**

**5 Digest DNA**

Add **95 µL rDNase** onto the column. Incubate at **room temperature** for **15 min**.



**+ 95 µL  
rDNase**

*Note: Centrifugation after incubation is not necessary.*

**RT, 15 min**

**6 Wash and dry silica membrane**

**1<sup>st</sup> wash**

Add **200 µL Buffer RB2** to the NucleoSpin® RNA Blood Column. Centrifuge for **30 s** at **11,000 x g**. Discard flow-through and Collection Tube and place the column into a new Collection Tube (2 mL; provided).



**+ 200 µL RB2**

*Buffer RB2 will inactivate the rDNase.*



**11,000 x g,  
30 s**

**2<sup>nd</sup> wash**

Add **600 µL Buffer RB3** to the NucleoSpin® RNA Blood Column. Centrifuge for **30 s** at **11,000 x g**. Discard flow-through and place the column into a new Collection Tube (2 mL; provided).



**+ 600 µL RB3**

*Note: Make sure that residual buffer from the previous steps is washed away with Buffer RB3, especially if the lysate has been in contact with the inner rim of the column during loading of the lysate onto the column. For efficient washing of the inner rim flush it with Buffer RB3.*



**11,000 x g,  
30 s**

**3<sup>rd</sup> wash**

Add **250 µL Buffer RB3** to the NucleoSpin® RNA Blood Column. Centrifuge for **2 min** at **11,000 x g**. In this step, ethanol is removed from the column.



**+ 250 µL RB3**

**11,000 x g,  
2 min**

Place the column into a nuclease-free Collection Tube (1.5 mL, supplied) and discard the Collection tube with flow-through from the previous step.

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

---

**7 Elute RNA**

Add **60 µL RNase-free H<sub>2</sub>O** (supplied) onto the column and centrifuge **30 s** at **11,000 x g**. The RNA is eluted into the Collection Tube.



**60 µL RNase-  
free H<sub>2</sub>O**

**11,000 x g,  
30 s**

For alternative elution procedures, see section 2.4.

---

## 5.2 RNA isolation from 400 µL blood

Before starting the preparation:

- Check if Wash Buffer RB3 and rDNase were prepared according to section 3.
- The complete procedure should be performed at room temperature (18–25 °C).
- The use of blood collected in common blood collection tubes with anticoagulant (typically EDTA) is recommended. For frozen blood samples, see section 2.3.
- Check if 70 % ethanol is available to adjust binding conditions.
- Refer to section 2.2 if less than 400 µL whole blood is used.

### 1 Lyse blood

Provide **400 µL whole blood** in a Collection Tube (2 mL, with lid; provided).



**400 µL blood**

Add **400 µL Lysis Buffer DL** to the tube and close the lid. Mix. If necessary, shortly spin to clean the lid.

**+ 400 µL DL**

Add **10 µL Liquid Proteinase K** and close the lid.

**+ 10 µL Liquid Proteinase K**

Incubate **3–15 min** at **room temperature** (18–25 °C) vigorously shaking the tube on a shaker (e.g., Eppendorf Thermoshaker, 1,400 rpm).

**RT, 3–15 min**

Centrifuge briefly to clean the lid (~ 1 s at ~ **2,000 x g**). Short spin only!



**~ 2,000 x g,  
~ 1 s**

### 2 Adjust RNA binding conditions

Add **400 µL 70 % ethanol** to the tube and mix vigorously.



**+ 400 µL  
70 % ethanol**

*Note: It is important to thoroughly mix the ethanol with the lysate. Recommended: Place tubes in a rack with lid. Close the rack lid and strongly shake the assembly. Alternatively, pipette the solution up and down ~ 5 times.*

**Mix**

Centrifuge briefly to clean the lid (~ 1 s at ~ **2,000 x g**). Short spin only!



**~ 2,000 x g,  
~ 1 s**



### 3 Bind RNA

Transfer **610 µL** and transfer lysate into a **NucleoSpin® RNA Blood Column** placed in a Collection Tube.



**Load 610 µL lysate**

*Note: Do not pipette more than 650 µL into the spin column, this will cause the column to overflow! Avoid formation of foam and aerosols! Avoid wetting the rim (edge) of the column.*

Centrifuge **30 s** at **11,000 x g**. Discard flow-through and Collection Tube. Place the column in a new Collection Tube (2 mL; provided).



**11,000 x g,  
30 s**

**Apply the remaining lysate** into the NucleoSpin® RNA Blood Column.



**Load residual lysate**

*Note: Do not pipette more than 650 µL into the spin column, this will cause the column to overflow! Avoid foam and aerosol formation! Avoid wetting the rim (edge) of the column.*

Centrifuge **30 s** at **11,000 x g**. Discard flow-through and Collection Tube. Place the column in a new Collection Tube (2 mL; provided).



**11,000 x g,  
30 s**

### 4 Desalt silica membrane

Add **350 µL MDB** (Membrane Desalting Buffer) onto the column and centrifuge **30 s** at **11,000 x g**.



**+ 350 µL MDB**

*Note: After centrifugation, the column can remain in the Collection Tube including the flow-through! The flow-through may be slightly brown. The flow-through can remain in the tube without disturbing DNA digestion.*



**11,000 x g,  
30 s**

### 5 Digest DNA

Add **95 µL rDNase** onto the column. Incubate at **room temperature** for **15 min**.



**+ 95 µL rDNase**

*Note: Centrifugation after incubation is not necessary.*

**RT, 15 min**

**6 Wash and dry silica membrane**

**1<sup>st</sup> wash**

Add **200 µL Buffer RB2** to the NucleoSpin® RNA Blood Column. Centrifuge for **30 s** at **11,000 x g**. Discard flow-through and Collection Tube and place the column into a new Collection Tube (2 mL; provided).

*Buffer RB2 will inactivate the rDNase.*



+ 200 µL RB2

**11,000 x g,  
30 s**

**2<sup>nd</sup> wash**

Add **600 µL Buffer RB3** to the NucleoSpin® RNA Blood Column. Centrifuge for **30 s** at **11,000 x g**. Discard flow-through and place the column into a new Collection Tube (2 mL; provided).

*Note: Make sure that residual buffer from the previous steps is washed away with Buffer RB3, especially if the lysate has been in contact with the inner rim of the column during loading of the lysate onto the column. For efficient washing of the inner rim flush it with Buffer RB3.*



+ 600 µL RB3

**11,000 x g,  
30 s**

**3<sup>rd</sup> wash**

Add **250 µL Buffer RB3** to the NucleoSpin® RNA Blood Column. Centrifuge for **2 min** at **11,000 x g**. In this step, ethanol is removed from the column.

Place the column into a nuclease-free Collection Tube (1.5 mL, supplied) and discard the Collection tube with flow-through from the previous step.

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



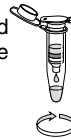
+ 250 µL RB3

**11,000 x g,  
2 min**

**7 Elute RNA**

Add **60 µL RNase-free H<sub>2</sub>O** (supplied) onto the column and centrifuge **30 s** at **11,000 x g**. The RNA is eluted into the Collection Tube.

For alternative elution procedures, see section 2.4.



**60 µL RNase-free  
H<sub>2</sub>O**

**11,000 x g,  
30 s**

## 6 NucleoSpin® RNA Blood Midi protocol – RNA isolation from 1.3 mL blood

Before starting the preparation:

- Check if Wash Buffer RB3 and rDNase were prepared according to section 3.
- The complete procedure should be performed at room temperature (18–25 °C).
- The use of blood collected in common blood collection tubes with anticoagulant (typically EDTA) is recommended. For frozen blood samples, see section 2.3.
- Check if 70 % ethanol is available to adjust binding conditions.
- For centrifugation, a centrifuge with a swing-out rotor and appropriate buckets capable of reaching 4,500 x g is required.
- Refer to section 2.2 if less than 1.3 mL whole blood is used.

### 1 Lyse blood

Provide **1.3 mL whole blood** in a 15 mL tube (provided).

Add **1.3 mL Lysis Buffer DL** to the tube and close the lid. Mix. If necessary, shortly spin to clean the lid.

Add **33 µL Liquid Proteinase K** and close the lid.

Incubate **3–15 min** at **room temperature** (18–25 °C) vigorously shaking the tube.

Centrifuge briefly to clean the lid (~ **1 s** at ~ **2,000 x g**). Short spin only!



**1.3 mL blood**  
**+ 1.3 mL DL**  
**+ 33 µL Liquid Proteinase K**  
**RT, 3–15 min**  
**~ 2,000 x g,**  
**~ 1 s**



### 2 Adjust RNA binding conditions

Add **1.3 mL 70 % ethanol** to the tube and mix vigorously.

*Note: It is important to thoroughly mix the ethanol into the lysate. Recommended: Vigorously shake for 5 s (e.g., on a vortexer at medium speed). Alternatively, pipette the solution up and down ~ 5 times.*

If necessary, centrifuge briefly to clean lid (~ **1 s** at ~ **2,000 x g**). Short spin only!

*This centrifugation step can be omitted if the lid is not wetted by the lysate. For example, mix by vortexing at medium speed or by pipetting up and down.*



**+ 1.3 mL**  
**70 % ethanol**  
**Mix**



**~ 2,000 x g,**  
**~ 1 s**

**3 Bind RNA**

Transfer the complete lysate (~ **4000 µL**) into a NucleoSpin® RNA Blood Midi Column placed in a 15 mL Collection Tube.

*Do not pipette more than 4000 µL into the Midi column, this will cause the column to overflow! Avoid foam and aerosol formation!*

Centrifuge **3 min** at **4,500 x g** and leave the column in the tube with the flow-through.



**Load max.  
4000 µL lysate**

**4,500 x g,  
3 min**

---

**4 Desalt silica membrane**

Add **1.2 mL MDB (Membrane Desalting Buffer)** onto the column and centrifuge **3 min** at **4,500 x g**.

Discard flow-through and Collection Tube and place the column in a new Collection Tube (15 mL, provided).



**+ 1.2 mL MDB**

**4,500 x g,  
3 min**

---

**5 Digest DNA**

Add **240 µL rDNase** onto the column.

Incubate at **room temperature** for **15 min** (centrifugation after this incubation is not necessary).



**+ 240 µL  
rDNase**

**RT,  
15 min**

---

**6 Wash and dry silica membrane****1<sup>st</sup> wash**

Add **1 mL Buffer RB2** to the NucleoSpin® RNA Blood Midi Column. Centrifuge for **3 min** at **4,500 x g**.

Leave the NucleoSpin® RNA Blood Midi Column in the tube with the flow-through.

**+ 1 mL RB2****4,500 x g,  
3 min****2<sup>nd</sup> wash**

Add **3 mL Buffer RB3** to the NucleoSpin® RNA Blood Midi Column. Centrifuge for **3 min** at **4,500 x g**.

Place the column into a nuclease-free Collection Tube (15 mL; provided) and discard the Collection Tube with flow-through from the previous step.

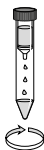
**+ 3 mL RB3****4,500 x g,  
3 min**

**Caution!** Ethanol carry-over may impair downstream analysis (e.g.,  $C_p$  retardation)! Therefore, removal of the column tube assembly from the centrifuge, transport to the lab bench, placement of it onto the bench and removal of the column from the tube should be done with great care in order to avoid any contamination of the column outlet with the flow-through/ethanol. To avoid any risk of wash buffer/ethanol contamination of the sample, discard wash buffer flow-through from the tube, re-insert the column into the tube and perform a further centrifugation (1–3 min 4,500 x g) in order to securely dry the column. Alternatively, use a fresh collection tube (not provided) for a secure column drying centrifugation.

**7 Elute RNA**

Add **200 µL RNase-free H<sub>2</sub>O** (supplied) onto the column. Centrifuge for **3 min** at **4,500 x g**. The RNA is eluted into the Collection Tube.

An additional elution step with **200 µL** fresh elution buffer will increase the total yield by approximately 25 %.

**+ 200 µL  
RNase-free  
H<sub>2</sub>O****4,500 x g,  
3 min**

## 7 Appendix

### 7.1 rDNase digestion in solution

The on-column rDNase digestion in the standard protocol is already very efficient and results in minimal residual DNA. This DNA will not be detectable in most downstream applications. Despite this, there are still certain applications which require even lower contents of residual DNA. However, removal of DNA to a completely undetectable level is challenging and the efficiency of an on-column DNA digestion is sometimes not sufficient for downstream applications requiring lowest residual content of DNA.

A typical example for such a demanding application is an RT-PCR reaction in which the primer molecules do not differentiate between cDNA (derived from RNA) and contaminating genomic DNA. Especially, if

- high copy number targets are analyzed (e.g., multi gene family, mitochondrial, plastidial 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 re-purification of the RNA (in order to remove buffer, salts, rDNase, and digested DNA) are usually required.

The high quality, recombinant, RNase-free DNase (rDNase) in the NucleoSpin® RNA Blood kits can also be used for a digestion in solution in order to remove trace amounts of contaminating DNA.

---

**1 Digest DNA (Reaction setup)**

Add **0.5 µL rDNase** per **10 µL eluted RNA** and mix moderately.

Centrifuge briefly (~ **1 s** at ~ **2,000 x g**) to collect all liquid in the lower part of the tube.

*Note: This step is important to ensure that every droplet of the RNA comes into contact with the rDNase to ensure efficient DNA digestion.*

---

**2 Incubate sample**

Incubate for **10 min** at **37 °C**.

---

**3 Repurify RNA**

Repurify RNA with a suitable RNA cleanup procedure, for example, using the NucleoSpin® RNA Clean-up kit (see ordering information) 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 H<sub>2</sub>O.

---

## 7.2 Troubleshooting

Problem	Possible cause and suggestions
RNA is degraded / no RNA obtained	<p><i>RNase contamination</i></p> <ul data-bbox="312 288 981 411" style="list-style-type: none"> <li>• Create an RNase-free working environment. Wear gloves during all steps of the procedure. Change gloves frequently. Use of sterile, disposable polypropylene tubes is recommended. Keep tubes closed whenever possible during the preparation. Glassware should be oven-baked for at least 2 hours at 250 °C before use.</li> </ul>
	<p><i>Reagents not applied or restored properly</i></p> <ul data-bbox="312 475 981 708" style="list-style-type: none"> <li>• Reagents not properly restored. Add the indicated volume of Reaction Buffer for rDNase and 96 % ethanol to Buffer RB3 Concentrate and mix. Reconstitute and store lyophilized rDNase according to instructions given in section 3.</li> <li>• Sample and reagents have not been mixed completely. Always vortex vigorously after each reagent has been added.</li> <li>• No ethanol has been added after lysis. Binding of RNA to the silica membrane is only effective in the presence of ethanol.</li> </ul>
Poor RNA quality or yield	<p><i>Kit storage</i></p> <ul data-bbox="312 767 953 948" style="list-style-type: none"> <li>• Reconstitute and store lyophilized rDNase according to instructions given in section 3.</li> <li>• Store other kit components at room temperature. Storage at low temperatures may cause salt precipitation.</li> <li>• Keep bottles tightly closed in order to prevent evaporation or contamination.</li> </ul>
	<p><i>Ionic strength and pH influence <math>A_{260}</math> absorption as well as ratio <math>A_{260}/A_{280}</math></i></p> <ul data-bbox="312 1007 981 1185" style="list-style-type: none"> <li>• For adsorption measurement, use 5 mM Tris pH 8.5 as diluent. Please see also: <ul data-bbox="351 1059 981 1185" style="list-style-type: none"> <li>- Manchester, K L. 1995. Value of <math>A_{260}/A_{280}</math> ratios for measurement of purity of nucleic acids. <i>Biotechniques</i> 19, 208–209.</li> <li>- Wilfinger, W W, Mackey, K and Chomczynski, P. 1997. Effect of pH and ionic strength on the spectrophotometric assessment of nucleic acid purity. <i>Biotechniques</i> 22, 474–481.</li> </ul> </li> </ul>



Problem	Possible cause and suggestions
Clogged NucleoSpin® Column / Poor RNA quality or yield	<i>Sample material</i>
	<ul style="list-style-type: none"> <li>• Bad sample quality. Make sure blood is collected into a standard blood collection tube (e.g., EDTA tube) according to the manufacturer's instructions; using fresh blood is always recommended. Sample should be stored at 4 °C for no longer than 24 hours. Freeze sample if it is not possible to process within one day.</li> </ul>
	<i>Inappropriate lysis / binding conditions</i>
	<ul style="list-style-type: none"> <li>• Do not premix Liquid Proteinase K with Lysis Buffer DL.</li> <li>• Make sure to vigorously shake during lysis incubation – shaking is essential for the procedure!</li> <li>• Make sure to use 70 % ethanol in this procedure to adjust binding conditions.</li> </ul>

Problem	Possible cause and suggestions
Contamination of RNA with genomic DNA	<p><i>rDNase not active</i></p> <ul style="list-style-type: none"> <li>• Reconstitute and store lyophilized rDNase according to instructions given in section 3.</li> </ul>
	<p><i>DNase solution not properly applied</i></p> <ul style="list-style-type: none"> <li>• Pipette rDNase solution directly onto the center of the silica membrane.</li> </ul>
	<p><i>High leukocyte number</i></p> <ul style="list-style-type: none"> <li>• The higher the leukocyte number, the higher the risk to detect residual DNA in the eluted RNA. To avoid this, use less blood.</li> </ul>
	<p><i>DNA detection system too sensitive</i></p> <ul style="list-style-type: none"> <li>• The amount of DNA contamination is effectively reduced during the on-column digestion with rDNase. However, it cannot be guaranteed that the purified RNA is 100 % free of DNA. Therefore, in very sensitive applications it might still be possible to detect DNA. The NucleoSpin® RNA system was checked by the following procedure: One million HeLa cells are subjected to RNA isolation. RNA eluate is used as a template for PCR detection of a 1 kb fragment in a 30 cycle reaction. Generally, no PCR product is obtained while skipping the DNase digest usually leads to positive PCR results.</li> <li>• The probability of DNA detection with PCR increases with: <ul style="list-style-type: none"> <li>- the number of DNA copies per preparation: single copy target &lt; plasmid / mitochondrial target &lt; plasmid transfected into cells</li> <li>- decreasing of PCR amplicon size.</li> </ul> </li> <li>• Use larger PCR targets (e.g., &gt; 500 bp) or intron spanning primers if possible.</li> </ul> <p><b>Use protocol 7.1 for subsequent rDNase digestion in solution.</b></p>
Suboptimal performance of RNA in downstream experiments	<p><i>Carry-over of ethanol or salt</i></p> <ul style="list-style-type: none"> <li>• Do not let the flow-through touch the column outlet after the second Buffer RB3 wash. Be sure to centrifuge at the corresponding speed for the respective time in order to remove ethanolic Buffer RB3 completely.</li> <li>• Check if Buffer RB3 has been equilibrated to room temperature before use. Washing at lower temperatures lowers efficiency of salt removal by Buffer RB3.</li> </ul>
	<p><i>Store isolated RNA properly</i></p> <ul style="list-style-type: none"> <li>• Eluted RNA should always be kept on ice for optimal stability since trace contaminations of omnipresent RNases (general lab ware, fingerprints, dust) will degrade the isolated RNA. For short term storage freeze at -20 °C, for long term storage freeze at -70 °C.</li> </ul>

### 7.3 Ordering information

Product	REF	Pack of
NucleoSpin® RNA Blood	740200.10 / .50	10 / 50
NucleoSpin® RNA Blood Midi	740210.20	20
NucleoSpin® 8 RNA Blood	740220 / .5	12 x 8 / 60 x 8
NucleoSpin® 96 RNA Blood	740225.2 / .4	2 x 96 / 4 x 96
NucleoSpin® miRNA Plasma	740981.10 / .50 / .250	10 / 50 / 250
NucleoSpin® RNA Clean-up	740948.10 / .50 / .250	10 / 50 / 250
NucleoSpin® RNA Clean-up XS	740903.10 / .50 / .250	10 / 50 / 250
NucleoSpin® RNA / DNA Buffer Set	740944	Suitable for 100 preps
NucleoSpin® RNA	740955.10 / .50 / .250	10 / 50 / 250
NucleoSpin® RNA Midi	740962.20	20
NucleoSpin® RNA / Protein	740933.10 / .50 / .250	10 / 50 / 250
NucleoSpin® TriPrep*	740966.10 / .50 / .250	10 / 50 / 250
NucleoSpin® miRNA	740971.10 / .50 / .250	10 / 50 / 250
NucleoSpin® RNA XS	740902.10 / .50 / .250	10 / 50 / 250
rDNase Set	740963	1 set
Collection Tubes (2 mL)	740600	1000

Visit [www.mn-net.com](http://www.mn-net.com) for more detailed product information.

### 7.4 Product use restriction/warranty

**NucleoSpin® RNA Blood** 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:**

NucleoSpin® is a trademark of MACHEREY-NAGEL GmbH & Co. KG

NucleoBond® is a trademark of MACHEREY-NAGEL GmbH & Co. KG.



[www.mn-net.com](http://www.mn-net.com)

## MACHEREY-NAGEL



MACHEREY-NAGEL GmbH & Co. KG · Neumann-Neander-Str. 6-8 · 52355 Düren · Germany

DE / International:

Tel.: +49 24 21 969-0  
 Fax: +49 24 21 969-199  
 E-mail: [info@mn-net.com](mailto:info@mn-net.com)

CH:

Tel.: +41 62 388 55 00  
 Fax: +41 62 388 55 05  
 E-mail: [sales-ch@mn-net.com](mailto:sales-ch@mn-net.com)

FR:

Tel.: +33 388 68 22 68  
 Fax: +33 388 51 76 88  
 E-mail: [sales-fr@mn-net.com](mailto:sales-fr@mn-net.com)

US:

Tel.: +1 484 821 0984  
 Fax: +1 484 821 1272  
 E-mail: [sales-us@mn-net.com](mailto:sales-us@mn-net.com)