



Genomic DNA from organs and cells

User manual

NucleoSpin[®] 96 DNA RapidLyse

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

For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Takara Bio Inc. ©2017

March 2017/Rev. 01

Contact MN

Germany and international

MACHEREY-NAGEL GmbH & Co. KG

Neumann-Neander-Str. 6–8 · 52355 Düren · Germany

Tel.: +49 24 21 969-0

Toll-free: 0800 26 16 000 (Germany only)

Fax: +49 24 21 969-199

E-mail: info@mn-net.com

Technical Support Bioanalysis

Tel.: +49 24 21 969-270

E-mail: tech-bio@mn-net.com

USA

MACHEREY-NAGEL Inc.

2850 Emrick Blvd. · Bethlehem, PA 18020 · USA

Tel.: +1 484 821 0984

Toll-free: 888 321 6224 (MACH)

Fax: +1 484 821 1272

E-mail: sales-us@mn-net.com

France

MACHEREY-NAGEL SARL à associé unique

1, rue Gutenberg · 67722 Hoerdt · France

Tel.: +33 388 68 22 68

Fax: +33 388 51 76 88

E-mail: sales-fr@mn-net.com

Switzerland

MACHEREY-NAGEL AG

Hirsackerstr. 7 · 4702 Oensingen · Switzerland

Tel.: +41 62 388 55 00

Fax: +41 62 388 55 05

E-mail: sales-ch@mn-net.com

www.mn-net.com

Table of contents

1 Components	4
1.1 Kit contents	4
1.2 Reagents to be supplied by user	4
1.3 About this user manual	4
2 Product description	5
2.1 The basic principle	5
2.2 Kit specifications	5
2.3 Required hardware	6
2.4 Automated processing on robotic platforms	8
2.5 Handling, preparation, and storage of starting materials	9
2.6 Lysis of sample material	9
2.7 Elution procedures	9
3 Storage conditions and preparation of working solutions	11
4 Safety instructions	12
5 Protocols	14
5.1 NucleoSpin® 96 DNA RapidLyse – centrifuge processing	14
5.2 NucleoSpin® 96 DNA RapidLyse – vacuum processing	19
5.3 Protocol for challenging samples (e.g., spleen and lung)	24
6 Appendix	26
6.1 Troubleshooting	26
6.2 Ordering information	28
6.3 Product use restriction/warranty	29

1 Components

1.1 Kit contents

REF	NucleoSpin® 96 DNA RapidLyse	
	1 x 96 preps 740110.1	4 x 96 preps 740110.4
Lysis Buffer RLY	30 mL	70 mL
Binding Buffer RLB	2 x 25 mL	2 x 100 mL
Wash Buffer RLW (Concentrate) ¹	50 mL	200 mL
Elution Buffer RLE ²	13 mL	60 mL
Liquid Proteinase K	1250 µL	4,5 mL
NucleoSpin® 96 DNA RapidLyse Binding Plates (green rings)	1	4
Square-well Blocks with self-adhering PE Foil	2	8
MN Wash Plates ³	1	4
User manual	1	1

1.2 Reagents to be supplied by user

- 96–100 % ethanol (for preparation of working solutions; see section 3)

For more detailed information regarding special hardware required for centrifuge or vacuum processing, please see section 2.3.

1.3 About this user manual

It is strongly recommended for first time users to read the detailed protocol sections of the **NucleoSpin® 96 DNA RapidLyse** kit before using this product. Experienced users, however, may refer to the Protocol-at-a-glance instead. 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 online at www.mn-net.com.

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

¹ For preparation of working solutions and storage conditions, see section 3.

² Elution Buffer RLE: 5 mM Tris/HCl, pH 8.5

³ For use with vacuum only.

2 Product description

2.1 The basic principle

The **NucleoSpin® 96 DNA RapidLyse** kit is designed for the efficient isolation of high molecular weight genomic DNA from cells and organs like liver, kidney, heart, muscle, spleen, and lung. Processing of mouse tail and ear clippings is also possible. Fresh, frozen, and ethanol-preserved samples can be used.

The **NucleoSpin® 96 DNA RapidLyse** kit lyses samples in maximal one hour agitated incubation at 56 °C. This is enabled by a thoroughly designed lysing setup with well balanced parameters that comprise a special lysis buffer in combination with Liquid Proteinase K. Incubation over night or for several hours is not necessary.

2.2 Kit specifications

Kit specifications at a glance

Parameter	NucleoSpin® 96 DNA RapidLyse
Technology	Silica-membrane technology
Format	96-well plates
Processing	Manual and automated, vacuum, positive pressure or centrifugation
Sample material	Fresh, frozen, dried, and ethanol preserved tissue samples (e.g., organs), eukaryotic cells
Sample amount	Up to 30 mg fresh weight (sample dependent)
Typical yield	Up to 4 µg DNA per mg tissue (sample dependent)
A_{260}/A_{280}	1.7–1.9
Elution volume	100 µL
Preparation time	60 min (96 preps, excluding lysis)
Lysis time	Maximal 1 h
Binding capacity	40 µg

- **NucleoSpin® 96 DNA RapidLyse** can be processed by vacuum, positive pressure or in a centrifuge and is designed for manual or automated use. The kit allows easy automation on common liquid handling instruments.
- The **NucleoSpin® 96 DNA RapidLyse** kit allows the purification of multiples of 96 samples. The kits are supplied with accessory plates for highest convenience. Further Accessory components (e.g., lysis plates, elution plates) can be individually selected from a variety of suitable accessories to provide highest flexibility (see section 2.3 for further information).

2.3 Required hardware

NucleoSpin® 96 DNA RapidLyse can be processed under vacuum, positive pressure or with centrifugation. Certain hardware for processing is required.

Centrifugation

For centrifugation, a microtiterplate centrifuge is required. This centrifuge must be able to accommodate the NucleoSpin® 96 DNA RapidLyse Binding Plate stacked on a Square-well Block and need to reach accelerations of 4,000 x *g* (bucket height: min. 75 mm).

Consumables for waste collection need to be ordered separately (e.g., MN Square-well Blocks). When processing two 96-well plates at once, six MN Square-well blocks are sufficient to collect the entire liquid waste without the need of emptying and reusing the MN Square-well Blocks (see ordering information, section 6.2). Alternatively, it is possible to use a single MN Square-well Block, which needs to be emptied after every centrifugation step.

Vacuum processing

The **NucleoSpin® 96 DNA RapidLyse** kit can be used with the NucleoVac 96 Vacuum Manifold. When using **NucleoSpin® 96 DNA RapidLyse** with less than 96 samples, Self-adhering PE Foil should be used in order to close and protect non-used wells of the NucleoSpin® 96 DNA RapidLyse Binding Plate and thus guarantee proper vacuum (see ordering information, section 6.2).

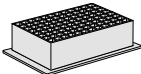
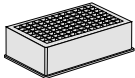
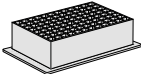
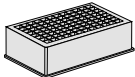

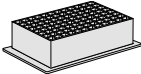
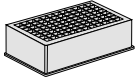
The manifold may be used with a vacuum pump, house vacuum, or water aspirator. We recommend a vacuum of -0.2 to -0.6 bar (reduction of atmospheric pressure). The use of the NucleoVac Vacuum Regulator (see ordering information, section 6.2) is recommended. Alternatively, adjust the vacuum in a way that the sample flows through the column with a rate of 1–2 drops per second. Depending on the amount of sample being used, the vacuum times may need to be increased for complete filtration.

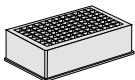
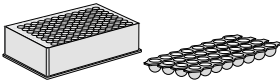

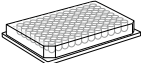
In addition, a suitable centrifuge for sample preparation steps may be required.

For general consumables and equipment needed, please see section 1.2.

Additional accessories for use of the NucleoSpin® 96 DNA RapidLyse Kit

Additional accessories for use of the **NucleoSpin® 96 DNA RapidLyse Kit** are available from MACHEREY-NAGEL (see ordering information, section 6.2).

Protocol step	Additional consumables	Remarks
Lyse samples	MN-Square-well Block	For sample lysis
		
	Square-well Block	
		
Adjust binding conditions	MN Square-well Block	Recommended for automated processing only
		
	Square-well Block	Self-adhering PE-Foil is required
		
Bind DNA	MN Wash Plate	MN Wash Plate minimizes the risk of cross contamination (vacuum processing only)
		
	MN Square-well Block	For waste collection during centrifugation (reusable)
		
	Square-well Block	
		

Protocol step	Additional consumables	Remarks
Elute DNA	Square-well Block 	
	Round-well Block with Cap Strips 	
	Rack of Tube Strips with Cap Strips 	
	Elution plates 	

2.4 Automated processing on robotic platforms

NucleoSpin® 96 DNA RapidLyse can be used fully automated on many common laboratory workstations. For the availability of scripts and general considerations about adapting the kit on a certain workstation, please contact MN. Full processing under vacuum or positive pressure enables complete automation without the need for centrifugation steps for drying of the binding membrane or for elution. However, a final elution step by centrifugation is recommended in order to achieve a DNA solution with higher concentration.

The risk of cross-contamination is reduced by optimized vacuum or positive pressure settings during the elution step and by the improved shape of the outlets of the NucleoSpin® 96 DNA RapidLyse Binding Plate.

Drying of the NucleoSpin® 96 DNA RapidLyse Binding Plates under vacuum or positive pressure is sufficient because the bottom of the plate is protected by the MN Wash Plate during the washing steps. Therefore it is recommended to integrate the MN Wash Plate into the automation procedure. The MN Frame (see ordering information, section 6.2) can be used to position the disposable MN Wash Plate inside the vacuum chamber. This also reduces the risk of cross-contamination, as common metal adaptors tend to get contaminated by gDNA. Thorough cleaning of the vacuum chamber is recommended after each run to prevent DNA-containing aerosols from forming.

Visit MN online at www.mn-net.com or contact your local MACHEREY-NAGEL distributor for technical support regarding hardware, software, setup instructions, and selection of the protocol.

2.5 Handling, preparation, and storage of starting materials

Fresh, frozen, and ethanol preserved samples can be used. Make sure not to use more than 30 mg sample.

2.6 Lysis of sample material

In order to obtain optimal DNA yields and a smooth processing, sample material should be thoroughly lysed. Most samples can be processed according to procedure 5.1. However, some sample materials (e.g., spleen or lung) need to be lysed according to an alternative protocol, which requires additional material (see section 5.2 and 6.2).

2.7 Elution procedures

It is possible to adjust the elution method and the volume of the elution buffer to the subsequent application of interest. In addition to the standard method described in the protocols (recovery rate about 70–90%), there are several modifications possible. Use elution buffer preheated at 70 °C for one of the following procedures:

- **High yield:** Perform two elution steps with the volume indicated in the individual protocol. About 90–100 % of bound nucleic acids can be eluted.
- **High concentration:** Perform one elution step with only 60 % of the volume indicated in the individual protocol. Concentration of DNA will be about 30 % higher than with the standard elution procedure. Maximum yield of bound nucleic acids is about 80 %.
- **High yield and high concentration:** Apply half the volume of elution buffer as indicated in the individual protocol, incubate for 3 min and centrifuge. Apply a second aliquot of elution buffer, incubate and centrifuge again. Thus, about 85–100 % of bound nucleic acids are eluted in the standard elution volume at a high concentration.
- **Convenient elution:** For convenience, elution buffer of ambient temperature may be used. This will result in a slightly lower yield (approximately 20 %) compared to elution with heated elution buffer.

Elution may also be performed with Tris-EDTA-buffer (TE) of pH equal or higher than 8. This will increase DNA stability during long term or multi-use storage at 4 °C (or ambient temperature) by inhibiting omnipresent DNases. However, high EDTA concentrations may interfere with certain downstream applications.

For optimal performance of isolated DNA in downstream applications, we recommend eluting with the supplied Elution Buffer RLE. For long term storage, the Elution Buffer RLE should be kept at - 20 °C. Several freeze-thaw cycles will not interfere with most downstream applications.

Performance of long-range PCR (e.g., > 10 kb) or the detection limit of trace amount of DNA species, may be reduced after multiple freeze-thaw cycles or prolonged storage of eluted DNA at 4 °C or room temperature. This is due to shearing of DNA or adsorption to surfaces.

Due to the dead volume of the silica membrane, please note that the difference between the dispensed elution buffer volume and the recovered elution buffer volume containing genomic DNA is approximately 40 μ L (recovered elution volume = dispensed elution volume - 40 μ L).

3 Storage conditions and preparation of working solutions

Attention: Buffer RLB contains chaotropic salts! Wear gloves and goggles!

CAUTION: Buffer RLB contains chaotropic 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.

Storage conditions:

- All kit components can be stored at room temperature (18–25 °C) and are stable for at least one year.
- Storage at lower temperatures may cause precipitation of salts. If a salt precipitation is observed, incubate the bottle at 30–40 °C for some minutes and mix well until all of the precipitate is redissolved. The performance of the kits is not affected by the salt precipitates.

Prior to the **NucleoSpin® 96 DNA RapidLyse** procedure, prepare the following:

- **Wash Buffer RLW:** Add the indicated volume of ethanol (96–100%) to Wash Buffer RLW Concentrate. Mark the label of the bottle to indicate that ethanol was added. **Wash Buffer RLW can be stored** at room temperature (18–25 °C) for up to one year.
- **Liquid Proteinase K** is ready to use. After first time use, store **Liquid Proteinase K** at 4 °C or -20 °C.

NucleoSpin® 96 DNA RapidLyse		
REF	1 x 96 preps 740110.1	4 x 96 preps 740110.4
Wash Buffer RLW (Concentrate)	50 mL Add 200 mL ethanol	200 mL Add 800 mL ethanol

4 Safety instructions

The following components of the **NucleoSpin® 96 DNA RapidLyse** 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>
RLB	Guanidinium thiocyanate 30–60 % and ethanol 5–20 % <i>Guanidinhydrochlorid 30–60 % und Ethanol 5–20 %</i> CAS 593-84-0, 64-17-5	 WARNING ACHTUNG	226, 302, 412, EUH031	210, 233, 260, 273, 301+312, 330, 370+378, 403+235
Proteinase K	Proteinase K, liquid 1–3 % <i>Proteinase K flüssig 1–3 %</i> CAS 39450-01-6	 WARNING ACHTUNG	317	261, 272, 280, 302+352, 333+313, 363

Hazard phrases

H226	Flammable liquid and vapour. <i>Flüssigkeit und Dampf entzündbar.</i>
H302	Harmful if swallowed. <i>Gesundheitsschädlich bei Verschlucken.</i>
H317	May cause an allergic skin reaction. <i>Kann allergische Hautreaktionen verursachen.</i>
H412	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

P210	Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking. <i>Von Hitze, heißen Oberflächen, Funken, offenen Flammen sowie anderen Zündquellenarten fernhalten. Nicht rauchen.</i>
P233	Keep container tightly closed. <i>Behälter dicht verschlossen halten.</i>
P260	Do not breathe dust/fume/gas/mist/vapours/spray. <i>Staub/Rauch/Gas/Nebel/Dampf/Aerosol nicht einatmen.</i>

- P261 Avoid breathing dust/fume/gas/mist/vapours/spray.
Einatmen von Staub/Rauch/Gas/Nebel/Dampf/Aerosol vermeiden.
- P272 Contaminated work clothing should not be allowed out of the workplace.
Kontaminierte Arbeitskleidung nicht außerhalb des Arbeitsplatzes tragen.
- P273 Avoid release to the environment.
Freisetzung in die Umwelt vermeiden.
- P280 Wear protective gloves/protective clothing/eye protection/face protection.
Schutzhandschuhe/Schutzkleidung/Augenschutz/Gesichtsschutz tragen.
- P301+312 IF SWALLOWED: Call a POISON CENTER/ doctor/.../ if you feel unwell.
BEI VERSCHLUCKEN: Bei Unwohlsein GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.
- P302+352 IF ON SKIN: Wash with plenty of water/...
BEI BERÜHRUNG MIT DER HAUT: Mit viel Wasser/... waschen.
- P330 Rinse mouth.
Mund ausspülen.
- P333+313 If skin irritation or rash occurs: Get medical advice/attention.
Bei Hautreizung oder -ausschlag: Ärztlichen Rat einholen/ärztliche Hilfe hinzuziehen.
- P363 Wash contaminated clothing before reuse.
Kontaminierte Kleidung vor erneutem Tragen waschen.
- P370+378 In case of fire: Use ... to extinguish.
Bei Brand: ... zum Löschen verwenden.
- P403+235 Store in a well-ventilated place. Keep cool.
An einem gut belüfteten Ort aufbewahren. Kühl halten.

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



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.

5 Protocols

5.1 NucleoSpin® 96 DNA RapidLyse – centrifuge processing

- For hardware requirements, refer to section 2.3.
- For storage conditions, refer to section 3.
- For use of recommended accessories, refer to section 2.4.

Before starting the preparation:

- Check if Buffer RLW was prepared according to section 3.
- Set incubator or oven to 56 °C.
- Set incubator or oven to 70 °C after lysis step.

1 Lyse sample

Place the samples into the wells of the **Square-well Block** or an appropriate lysis container (e.g. U-bottom plate or 2 mL tube).

Add **150 µL Buffer RLY** to each sample.

+ 150 µL RLY

Add **10 µL Liquid Proteinase K** to each sample.

+ 10 µL Liquid Proteinase K

Seal the plate with a Self-adhering Foil and **mix** by vigorous shaking for 10–15 s. Spin briefly (15 s; 1,500 x g) to collect any sample at the bottom of the wells.

Mix

Incubate at **56 °C** on a thermomixer at maximum speed until the sample appears visually lysed.

**56 °C,
max. 1 h**

Note: Lysis time depends on sample material and may vary from a couple of minutes up to one hour (for mammalian cells reduce incubation to 30 min).

Do not use conical plates or 1.5 mL tubes, as their shape will impair thorough mixing. Use the Square-well Block or 2 mL tubes which will facilitate proper sample and lysis buffer agitation. Make sure that the tissue sample is submerged in the lysis buffer!

Take care not to moisten the rims of the individual wells while dispensing the buffer.

Spin briefly the Square-well Block or lysis container (15 s; 1,500 x g) to collect any condensate at the bottom of the wells.

Residual hair and/or bones in the lysate can be removed by centrifugation (2 min; 5,600–6,000 x g) and transfer of the supernatant to a new Square-well Block (not supplied with the kit).

2 Adjust DNA binding conditions

Add **440 µL Buffer RLB to each sample** and **mix** (e.g., pipetting up and down, pulse vortexing or shaking). Seal the plate with a Self-adhering Foil if necessary.

**+ 440 µL RLB
Mix**

Take care not to moisten the rims of the individual wells while dispensing the buffer. Spin briefly the Square-well Block or lysis container (15 s; 1,500 x g) to collect any condensate at the bottom of the wells.

Never centrifuge at higher g-forces or for longer periods as DNA will precipitate.

Place a NucleoSpin® 96 DNA RapidLyse Binding Plate on a Square-well Block.

If using more than one plate, label the plates for later identification. The use of a second plate placed on a MN Square-well Block avoids the need to balance the centrifuge.

3 Transfer lysates

Transfer the **lysates** resulting from step 3 carefully from the Square-well Block into the wells of the **NucleoSpin® 96 DNA RapidLyse Binding Plate**.

Transfer lysate

Take care to dispense the lysate centric onto the silica membrane.

Do not moisten the rims of the individual wells while dispensing the samples – moistened rims may cause cross contamination during centrifugation.

We recommend the use of an electronic eight-channel pipetting device with extra long tips capable of holding approx. 1000 µL, for transfer of the lysate from the Square-well Block to the NucleoSpin® 96 DNA RapidLyse Binding Plate.

4 Bind DNA

After transfer, seal the openings of the plate with Self-adhering PE-Foil and incubate the lysates for **1 min** on the NucleoSpin® 96 DNA RapidLyse Binding Plate at **room-temperature**.

1 min, RT

Place the Square-well Blocks with NucleoSpin® 96 DNA RapidLyse Binding Plates onto the centrifuge carriers and insert them into the rotor buckets. Centrifuge for **2 min** at **4,000 x g**.

**4,000 x g,
2 min**

Typically, the lysates will have passed through the silica membrane within a few minutes. The centrifugation process can be extended to 20 min, if the lysates have not passed completely.

Discard Square-well block with flow through. Put NucleoSpin® 96 DNA RapidLyse Binding Plate onto a new Square-well Block (not provided).

When using NucleoSpin® 96 DNA RapidLyse with less than 96 samples, Self-adhering PE Foil should be used in order to close and protect non-used wells of the NucleoSpin® 96 DNA RapidLyse Binding Plate and thus guarantee proper vacuum.

5 Wash silica membrane**1st wash**

Remove the Self-adhering PE Foil and add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Seal the plate with a new Self-adhering PE Foil and centrifuge again for **2 min** at **4,000 x g**.

+ 750 µL RLW
4,000 x g,
2 min

Discard flow through and place NucleoSpin® 96 DNA RapidLyse Binding Plate back onto the Square-well Block.

2nd wash

Remove the Self-adhering PE Foil and add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Seal the plate with a new Self-adhering PE Foil and centrifuge again for **2 min** at **4,000 x g**.

+ 750 µL RLW
4,000 x g,
2 min

Discard flow through and place NucleoSpin® 96 DNA RapidLyse Binding Plate back onto the Square-well Block.

3rd wash

Remove the Self-adhering PE Foil and add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Seal the plate with a new Self-adhering PE Foil and centrifuge again for **2 min** at **4,000 x g**.

+ 750 µL RLW
4,000 x g,
2 min

Discard flow through and place NucleoSpin® 96 DNA RapidLyse Binding Plate back onto the Square-well Block (not provided).

6 Dry silica membrane

Remove the Self-adhering PE Foil and place the NucleoSpin® 96 DNA RapidLyse Binding Plate onto a Square-well Block. Place it in an incubator for **10 min** at **70 °C** to evaporate residual ethanol.

**70 °C,
10 min**

Removal of ethanol by evaporation at 70 °C is more effective than prolonged centrifugation.

or

Optional:

Centrifuge for **5 min** at **4,000 x g**.

**4,000 x g,
5 min**

Note: Residual wash buffer is removed in this step.

7 Elute highly pure DNA

Place the NucleoSpin® 96 DNA RapidLyse Binding Plate onto a fresh Square-well Block and dispense **100 µL Buffer RLE** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Dispense the buffer directly onto the membrane. Incubate at room temperature for **1 min**.

+ 100 µL RLE

1 min, RT

Centrifuge for **5 min** at **4000 x g**.

If elution in small volume tubes is desired, place a 96 PCR plate (not supplied) on top of a Round-well Block or a Rack of Tube Strips and elute into the PCR plate.

**4,000 x g,
5 min**

Note: DNA yield can be increased by incubation for 4 min at room temperature before centrifugation.

For alternative elution procedures see section 2.5.

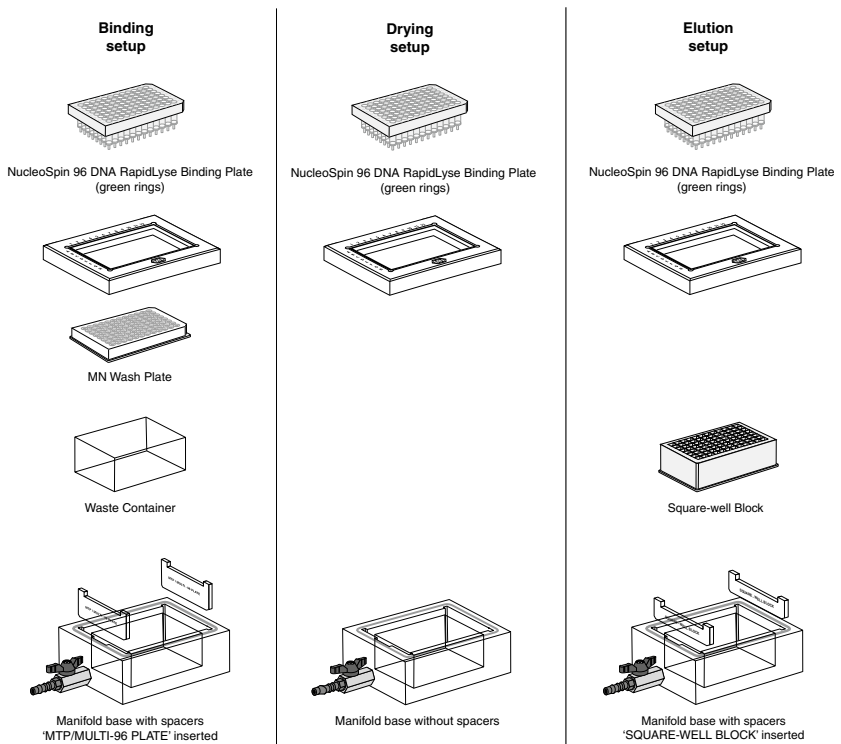
Setup of vacuum manifold:

5.2 NucleoSpin® 96 DNA RapidLyse – vacuum processing

- For hardware requirements, refer to section 2.3.
- For storage conditions, refer to section 3..
- For use of recommended accessories, refer to section 2.4.

Before starting the preparation:

- Check if Buffer RLW was prepared according to section 3.
- Set incubator or oven to 56 °C.



1 Lyse sample

Place the samples into the wells of the **Square-well Block** or an appropriate lysis container (e.g. U-bottom plate or 2 mL tube).

Add **150 µL Buffer RLY** to each sample.

+ 150 µL RLY

Add **10 µL Liquid Proteinase K** to each sample.

+ 10 µL Liquid Proteinase K

Seal the plate with a Self-adhering Foil and **mix** by vigorous shaking for 10–15 s. Spin briefly (15 s; 1,500 x g) to collect any sample at the bottom of the wells.

Mix

Incubate at **56 °C** on a thermomixer at maximum speed until the sample appears visually lysed.

Note: Lysis time depends on sample material (for mammalian cells reduce incubation to 30 min) and may vary from a couple of minutes up to one hour.

**56 °C,
max. 1 h**

Do not use conical plates or 1.5 mL tubes as their shape will impair thorough mixing. Use the Square-well Block or 2 mL tubes which will facilitate proper sample and lysis buffer agitation. Make sure that the tissue sample is submerged in the lysis buffer!

Take care not to moisten the rims of the individual wells while dispensing the buffer.

Spin briefly the Square-well Block or lysis container (15 s; 1,500 x g) to collect any condensate at the bottom of the wells.

Residual hair and/or bones in the lysate can be removed by centrifugation (2 min; 5,600–6,000 x g) and transfer of the supernatant to a new Square-well block (not supplied with the kit).

2 Adjust DNA binding conditions

Add **440 µL Buffer RLB** to each sample and **mix** (e.g., pipetting up and down, pulse vortexing or shaking). Seal the plate with a Self-adhering Foil if necessary.

+ 440 µL RLB

Mix

Take care not to moisten the rims of the individual wells while dispensing the buffer. Spin briefly the Square-well Block or lysis container (15 s; 1,500 x g) to collect any condensate at the bottom of the wells.

3 Transfer lysates

Transfer the lysates resulting from step 3 carefully from the Square-well block into the wells of the **NucleoSpin® 96 DNA RapidLyse Binding Plate**. **Transfer lysates**

Take care to dispense the lysate centric onto the silica membrane.

Do not moisten the rims of the individual wells while dispensing the samples – moistened rims may cause cross contamination during centrifugation.

We recommend the use of an electronic eight-channel pipetting device with extra long tips capable of holding approx. 1000 µL, for transfer of the lysate from the Square-well block to the NucleoSpin® 96 DNA RapidLyse Binding Plate.

Seal non-used wells of the NucleoSpin® 96 DNA RapidLyse Binding Plate with a Self-adhering PE-foil

Prepare the NucleoVac 96 Vacuum Manifold:

Place waste tray into vacuum manifold base. Insert spacers labeled “MTP/Multi-96 plate” notched side up and place the MN Wash Plate on them. Close the manifold with the manifold lid and place a NucleoSpin® 96 DNA RapidLyse Binding Plate on top of the manifold.

4 Bind DNA

Incubate the lysates for 1 min on the **NucleoSpin® 96 DNA RapidLyse Binding Plate**. **1 min, RT**

Apply vacuum until all lysates have passed through the wells of the NucleoSpin® DNA RapidLyse Binding Plate **(-0.4 bar*, 2 min)**. Release the vacuum. **-0.4 bar*, 2 min**

* reduction of atmospheric pressure

5 Wash silica membrane**1st wash**

Add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Apply vacuum until all buffer has passed through the wells (**-0.4 bar***; **2 min**).
Release the vacuum.

+ 750 µL RLW
-0.4 bar*,
2 min

2nd wash

Add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Apply vacuum until all buffer has passed through the wells (**-0.4 bar***; **2 min**).
Release the vacuum.

+ 750 µL RLW
-0.4 bar*,
2 min

3rd wash

Add **750 µL Buffer RLW** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Apply vacuum until all buffer has passed through the wells (**-0.4 bar***; **2 min**).
Release the vacuum.

+ 750 µL RLW
-0.4 bar*,
2 min

Remove MN Wash Plate

After the final washing step, close the valve, release the vacuum and remove the NucleoSpin® 96 DNA RapidLyse Binding Plate. Put it on a clean paper towel to remove residual EtOH-containing wash buffer. Remove manifold lid, MN Wash Plate, and waste container from the vacuum manifold.

6 Dry silica membrane

Insert the NucleoSpin® 96 DNA RapidLyse Binding Plate onto the lid, and close the manifold (without waste container). Apply maximum vacuum (**at least -0.6 bar***) for **10 min** to dry the membrane completely.

-0.6 bar*,
10 min

Note: Residual wash buffer is removed in this step.

Finally, release the vacuum.

* reduction of atmospheric pressure

7 Elute highly pure DNA

Insert spacers “Square-well Block” into the NucleoVac Vacuum Manifold’s short sides. Place a Square-well Block onto the spacer. Close the vacuum manifold and place the NucleoSpin® 96 DNA RapidLyse Binding Plate on top. Dispense **100 µL Buffer RLE** to each well of the NucleoSpin® 96 DNA RapidLyse Binding Plate. Dispense the buffer directly onto the membrane. Incubate at room temperature for **1 min**.

+ 100 µL RLE

1 min, RT

Apply vacuum for elution (**-0.6 bar***; **5 min**). Release the vacuum and seal the Square-well block with Self-adhering Foil.

**-0.6 bar*,
5 min**

Spin the Square-well block briefly to collect all samples at the bottom of the plate.

For alternative elution procedures see section 2.5.

* reduction of atmospheric pressure

5.3 Protocol for challenging samples (e.g., spleen and lung)

- For hardware requirements, refer to section 2.3.
- For storage conditions, refer to section 3.
- For use of recommended accessories, refer to section 2.4.

Before starting the preparation:

- Check if Buffer RLW was prepared according to section 3.
- Set incubator or oven to 56 °C.

The following items are additionally required for this protocol:

- MN Bead Tube Holder, NucleoSpin® Bead Tubes Type F (see ordering information, section 6.2).

1 Lyse sample

Place the sample into a **NucleoSpin® Bead Tube Type F**.

Add **120 µL Buffer RLE**. **+ 120 µL RLE**

Add **20 µL Buffer RLB**. **+ 20 µL RLB**

Add **10 µL Proteinase K**. **+ 10 µL Liquid Proteinase K**

Insert the Bead Tube into the **MN Bead Tube Holder** and **shake 20 min** at **full speed** on a Vortex-Genie® 2. Up to 30 mg of wet weight sample can be processed.

Shake 20 min, full speed

Note: The use of other disruption devices is not recommended in conjunction with Bead Tube Type F. Due to the lysing matrix (corundum and steel balls) high impact disruption devices will cause steel abrasion and possible demolition of the bead tubes!

2 Adjust DNA binding conditions

Add **420 µL Buffer RLB** and **mix** (e.g., vortex 3 s). **+ 420 µL RLB**

Centrifuge the tube at **11,000 x g** for approx. **5 s** (short spin), in order to clean the lid and sediment the lysing matrix.

Mix
11,000 x g,
5 s

DO NOT centrifuge for longer times and/or higher g-force, as this might damage the Bead Tubes due to the high density of the steel balls.

3 Bind DNA

Load cleared **supernatant** (approximately 500 µL) on the wells of the **NucleoSpin® 96 DNA RapidLyse Binding Plate**.

Load supernatant

Note: Do not disturb the lysing matrix. Make sure not to transfer corundum matter from the lysing tube onto the column!

Take care to dispense the lysate centric onto the silica membrane.

Do not moisten the rims of the individual wells while dispensing the samples – moistened rims may cause cross contamination during centrifugation.

We recommend the use of an electronic eight-channel pipetting device with extra long tips capable of holding approx. 1000 µL, for transfer of the lysate from the Square-well block to the NucleoSpin® 96 DNA RapidLyse Binding Plate.

Follow step 4 (see section 5.1) of the vacuum processing protocol, starting with the preparation of the NucleoVac 96 Vacuum Manifold or step 4 (see section 5.2) of the centrifuge processing protocol starting.

6 Appendix

6.1 Troubleshooting

Problem	Possible cause and suggestions
No or poor DNA yield	<i>Incomplete lysis</i>
	Sample has not completely been submerged during heat incubation. Cut samples into small pieces. Mix well. Be sure that the samples are fully submerged in Buffer RLY / Proteinase K mixture.
	<i>Reagents not applied properly</i>
	Prepare Wash Buffer RLW according to instructions (section 3).
No or poor DNA yield	<i>Suboptimal elution of DNA from the column</i>
	Preheat Buffer RLE to 70 °C before elution. Apply Buffer RLE directly onto the center of the silica membrane.
	Elution efficiencies decrease dramatically if elution is done with buffers with pH < 7. Use slightly alkaline elution buffer like Buffer RLE (pH 8.5).
	Perform a second elution step using 100 µL RLE and incubate for 5 min at RT before applying vacuum.
	<i>For alternative elution procedures see section 2.5.</i>
Poor DNA quality	<i>High A_{260}/A_{280} ratio</i>
	Ratios > 1.9 can be caused by RNA contamination. Usually, such RNA contamination does not interfere with downstream applications. Depending on sample type, amount, and disruption procedure, preparations might contain small amounts of RNA. If it is necessary to reduce RNA contamination to the lowest possible level, incubate the lysate after disruption for 5 min at 70 °C in order to inactivate the Proteinase K. After cooling to room temperature, add 20 µL RNase A (20 mg/mL) and incubate 5 min. Continue with the application of the lysate onto the silica membrane.
	<i>Reagents not applied properly</i>
	Prepare Buffer RLW according to instructions (see section 3).
	<i>Low A_{260}/A_{230} ratio</i>
	Increase the volume of Wash Buffer RLW to 900 µL per washing steps to remove residual impurities.

Problem	Possible cause and suggestions
RNA contamination	<p><i>RNA in sample</i></p> <p>If DNA free of RNA is desired, cool down to room temperature after lysis incubation and add 20 µL of an RNase A solution (20 mg/mL; see ordering information 6.2). Incubate for 15 min with moderate shaking.</p>
Poor performance of genomic DNA in enzymatic reactions	<p><i>Carry-over of ethanol</i></p> <p>After washing with Buffer RLW, centrifuge ≥ 4 min at 4000 x g in order to remove ethanolic Buffer RLW completely and evaporate residual ethanol by incubating the NucleoSpin® 96 DNA RapidLyse Binding Plate at 70 °C for 10 min.</p> <p>Increase vacuum drying time to 15 min.</p> <p><i>Contamination of DNA with inhibitory substances</i></p> <p>Do not elute DNA with TE buffer. EDTA may inhibit enzymatic reactions. Repurify DNA and elute in Buffer RLE.</p>
Clogged wells	<p><i>Too much sample material used</i></p> <p>Reduce the sample amount or follow procedure 5.2 for the next preparation.</p> <p><i>Hair or bones left in the lysate after step 2</i></p> <p>Centrifuge the Square-well Block for 3 min at 5,600–6,000 x g. Transfer lysates to a new Square-well Block without disturbing the debris pellet.</p> <p><i>Incomplete passage of lysate in step 4</i></p> <p>If no more than 300–500 µL of lysate is remaining in the columns, continue with step 5. Through the addition of Buffer RLW the sample is diluted and thus the sample will pass the column more easily.</p> <p><i>Poor vacuum pressure</i></p> <p>Check the vacuum manifold for a leaking air-flow. Press the NucleoSpin® DNA Binding Plate slightly onto the vacuum manifold lid before applying vacuum.</p>
Grayish lysate or membrane	<p><i>Lysis with Bead Tube Type F for 20 min on the MN Bead Tube Holder might cause a slight grayish color of the lysate, which is tolerable. Prolonged shaking or use of other disruption devices can cause steel abrasion.</i></p> <p>Do not perform prolonged incubation, do not use other disruption devices with Bead Tube Type F.</p>

6.2 Ordering information

Product	REF	Pack of
NucleoSpin® 96 DNA RapidLyse	740110.1	1 x 96 preps
	740110.4	4 x 96 preps
Liquid Proteinase K	740396	5 mL
RNase A	740505	50 mg
	740505.50	100 mg
Square-well Block	740481	4
	740481.24	24
Rack of Tube Strips (1 set consists of 1 rack, 12 strips with 8 tubes each, and 12 Cap Strips)	740477	4 sets
	740477.24	24 sets
Round-well Block with Cap Strips (1 set consists of 1 Round-well Block and 12 Cap Strips)	740475	4 sets
	740475.24	24 sets
MN Wash Plate	740479	4
	740479.24	24
Cap Strips	740478	48
	740478.24	288
NucleoVac 96 Vacuum Manifold	740681	1
NucleoVac Vacuum Regulator	740641	1
Self-adhering PE Foil	740676	50
Rubber Pad	740640	2
MN Bead Tube Holder	740469	1 piece
NucleoSpin® Bead Tubes Type F (1–3 mm corundum and 3 mm steel balls, recommended for challenging samples in combination with NucleoSpin® DNA RapidLyse – use only with MN Bead Tube Holder)	740816.50	50 pieces

Visit www.mn-net.com for more detailed product information.

6.3 Product use restriction/warranty

NucleoSpin® 96 DNA RapidLyse 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.

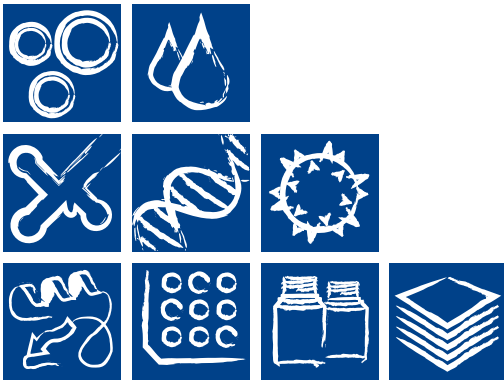
Product claims are subject to change. Therefore please contact our Technical Service Team for the most up-to-date information on MACHEREY-NAGEL products. You may also contact your local distributor for general scientific information. Applications mentioned in MACHEREY-NAGEL literature are provided for 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 (0) 24 21 969 270
e-mail: tech-bio@mn-net.com



Plasmid DNA
Clean-up
RNA
Genomic DNA
Viral RNA and DNA
Protein
High throughput
Accessories
Auxiliary tools



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

CH:

Tel.: +41 62 388 55 00
 Fax: +41 62 388 55 05
 E-mail: 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

US:

Tel.: +1 484 821 0984
 Fax: +1 484 821 1272
 E-mail: sales-us@mn-net.com