



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



takarabio.com

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

Genomic DNA from FFPE samples

User manual

NucleoSpin® 96 DNA FFPE

May 2015 / Rev. 01

MACHEREY-NAGEL

www.mn-net.com



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 EURL

1, rue Gutenberg · 67722 Hoerd · 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	Material to be supplied by user	4
2	Product description	5
2.1	The basic principle	5
2.2	Kit specifications	5
2.3	Handling, preparation, and storage of starting materials	6
2.4	Required hardware	7
2.5	Automated processing on robotic platforms	7
2.6	Elution procedures	8
3	Storage conditions and preparation of working solutions	9
4	Safety instructions	10
5	Protocols	12
5.1	Centrifuge processing	12
5.2	Vacuum processing	17
6	Appendix	23
6.1	Troubleshooting	23
6.2	Ordering information	25
6.3	Product use restriction/warranty	26

1 Components

1.1 Kit contents

NucleoSpin® 96 DNA FFPE		
REF	1 x 96 preps 740240.1	4 x 96 preps 740240.4
Lysis Buffer FL	30 mL	4 x 30 mL
Binding Buffer BQ1	45 mL	3 x 45 mL
Wash Buffer BW	100 mL	300 mL
Wash Buffer B5	100 mL	2 x 100 mL
Elution Buffer BE	30 mL	125 mL
Proteinase Buffer PB	15 mL	35 mL
Paraffin Dissolver (blue)	60 mL	2 x 125 mL
Decrosslink Buffer D-Link	30 mL	2 x 30 mL
Proteinase K* (lyophilized)	75 mg	4 x 75 mg
NucleoSpin® DNA FFPE Binding Plate	1	4
MN Wash Plate	1	4
Round-well Block (including Cap Strips)	1	4
MN Square-well Block	1	4
Rack of Tube Strips (including Cap Strips)	1	4
Self adhering PE Foil	5	10
User manual	1	1

1.2 Material to be supplied by user

Reagents

- 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 2.4.

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

2 Product description

Formalin-fixed, paraffin-embedded (FFPE) tissue samples are routinely prepared from human surgical tissue samples by fixation with formalin and embedding in paraffin. Thin sections of FFPE samples are commonly subjected to histopathological analysis and remaining paraffin-tissue blocks are usually archived. Existing extensive archives of FFPE tissue samples represent a valuable source for retrospective studies of gene expression patterns and mutation analysis. However, the use of such samples for DNA analysis is limited due to chemical modification by formaldehyde and fragmentation of the DNA during tissue processing (sampling, fixing, embedding) and storage (humidity, time, temperature) of the samples. Standard DNA isolation procedures often result in low DNA yield or poor performance in downstream applications (e.g., PCR).

2.1 The basic principle

The **NucleoSpin® 96 DNA FFPE** kit is designed for the isolation of DNA from formalin-fixed, paraffin-embedded (FFPE) tissue specimen. The procedure replaces the use of flammable and malodorous xylene or d-limonene commonly used for deparaffinization. Further, the procedure does not require the difficult removal of organic solvent from often barely visible tissue pellets. **NucleoSpin® 96 DNA FFPE** employs the odorless Paraffin Dissolver (patented) and allows effective lysis in a convenient two-phase system. This kit provides reagents for a convenient, reliable, and fast method to isolate DNA from 96 or 384 samples. First, the paraffin of FFPE sections is dissolved in the Paraffin Dissolver. Tissue is then digested by Proteinase K to solubilize the fixed tissue and release DNA into solution.

Appropriate conditions for binding of DNA to the silica membrane in the NucleoSpin® DNA FFPE Binding Plate are created by addition of large amounts of chaotropic salt and ethanol to the lysate. The binding process is reversible and specific to nucleic acids. While DNA is kept on the silica membrane, contaminations are removed by washing with two different wash buffers. Pure genomic DNA is finally eluted under low ionic strength conditions in a slightly alkaline elution buffer.

2.2 Kit specifications

The **NucleoSpin® 96 DNA FFPE** kit is recommended for the isolation of DNA from formalin-fixed, paraffin-embedded (FFPE) tissue samples. Samples are typically thin sections (approx. 3–20 µm thickness) of human or animal origin usually obtained by tissue resection or biopsy.

- **Sample amount:** The maximum sample size is determined by a) the amount of tissue and b) by the amount of paraffin. **NucleoSpin® 96 DNA FFPE** is suitable for up to 10 mg tissue. The amount of paraffin is limited to 15 mg (approx 17 µL), when using the standard protocol with Paraffin Dissolver. However, larger amounts of paraffin samples may be processed by using either additional Paraffin Dissolver or by deparaffinization using xylene. Typically, one or two sections of 10 µm to 20 µm thickness are sufficient for DNA purification. DNA yield depends on the amount or size of the embedded tissue samples. Excess of paraffin may be removed by trimming the paraffin sections or the paraffin embedded sample blocks before slicing.

- **DNA yield** strongly depends on the sample type, quality, quantity as well as conditions and duration of storage. Further, measured DNA yield may vary considerably between different quantification methods. Yield determined by absorption measurement at 260 nm or by a fluorescent dye (e.g., PicoGreen®) may deviate from values obtained by quantification with PCR. Even quantification values obtained via PCR with a short (e.g., 80 bp) and a long (e.g., 300 bp) amplicon may also differ considerably. The deviation of quantification also depends on DNA size distribution as well as on efficiency of decrosslinking (or extent of remaining crosslinks).
- **DNA size distribution:** DNA isolated from formalin-fixed, paraffin-embedded tissue shows a size distribution from 50 to 5,000 bases. Predominantly DNA of approx. 100–300 bases is observed, especially when the sample material is old. However, samples which were subjected to good tissue fixation, embedding, and storage conditions can yield DNA even larger than 5,000 bp.
- **DNA preparation time** strongly depends on the sample and the required lysis time. For best results lysis is performed at room temperature for at least three hours. For some kinds of samples a longer lysis (e.g., overnight) will even result in remarkably higher DNA yield.
- **NucleoSpin® 96 DNA FFPE** can be processed by vacuum or by centrifugation. The kit allows easy automation on common liquid handling instruments.
- The **NucleoSpin® 96 DNA FFPE** kits allow for the purification of multiples of 96 samples. The kits are supplied with accessory plates for highest convenience. The kits are designed for manual or automated use in a centrifuge or for use with a vacuum manifold.

2.3 Handling, preparation, and storage of starting materials

Many factors influence the yield and quality of DNA obtained from FFPE samples. The procedure of tissue sampling, post sampling delay before fixation, fixation time, embedding, and storage conditions have a high impact on DNA quality and yield. Starting from a paraffin embedded tissue block, samples should be sectioned under clean conditions. Paraffin sections may be stored at +4 °C or lower for at least several weeks without observable effects on DNA yield or usability. Long term storage of paraffin sections may have a negative effect on the DNA due to air oxidation. Wear gloves at all times during the preparation. Change gloves frequently.

2.4 Required hardware

NucleoSpin® 96 DNA FFPE can be processed under vacuum or by centrifugation. Certain hardware for processing is required.

Centrifugation

For centrifugation, a microtiterplate centrifuge is required. This centrifuge must be able to accommodate the NucleoSpin® DNA FFPE Binding Plate stacked on a Round- or Square-well Block and reach accelerations of 5,600–6,000 x *g* (bucket height: 85 mm). Regarding waste collection, suitable consumables (e.g., MN Square-well Blocks) are necessary and they are not included in the kit. For the most convenient handling, without the need of emptying and reusing the MN Square-well Blocks, it is recommended using six MN Square-well Blocks if two 96-well plates are processed at once (see ordering information, section 6.2). Alternatively, it is possible to empty the MN Square-well Blocks after every centrifugation step, reducing the amount of MN Square-well Blocks needed.

Vacuum processing

The **NucleoSpin® 96 DNA FFPE** kit can be used with the NucleoVac 96 Vacuum Manifold (see ordering information, section 6.2). When using **NucleoSpin® 96 DNA FFPE** with less than 96 samples, Self-adhering PE Foil (see ordering information, section 6.2) should be used in order to close and protect non-used wells of the NucleoSpin® DNA FFPE Binding Plate and thus guarantee proper vacuum. Establish a reliable vacuum source for the NucleoVac 96 Vacuum Manifold. The manifold may be used with a vacuum pump, house vacuum, or water aspirator. It is recommended a vacuum of -0.2 to -0.4 bar (reduction of atmospheric pressure). The use of the NucleoVac Vacuum Regulator (see ordering information, section 6.2) is recommended. Alternatively, adjust the vacuum so that during the purification 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. Additionally, a suitable centrifuge for sample preparation steps may be required.

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

2.5 Automated processing on robotic platforms

NucleoSpin® 96 DNA FFPE can be used fully automated on many common laboratory workstations. For the availability of scripts and general considerations about adapting **NucleoSpin® 96 DNA FFPE** on a certain workstation, please contact MN. Full processing under vacuum 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 higher concentrated eluted DNA. The risk of cross-contamination is reduced by optimized vacuum settings during the elution step and by the improved shape of the outlets of the NucleoSpin® DNA FFPE Binding Plate. Drying of the NucleoSpin® DNA FFPE Binding Plates under vacuum is sufficient because the bottom of the plate is protected by the MN Wash Plate during the washing steps. It is recommended to integrate the MN Wash Plate into the automated procedure to avoid cross-contamination by residual wash buffer. 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.

2.6 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, EDTA interferes, depending on the final concentration, with certain downstream applications.

For optimal performance of isolated DNA in downstream applications, it is recommended eluting with the supplied elution buffer and storing it, especially for long term 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 20 µL (recovered elution volume = dispensed elution volume - 20 µL).

3 Storage conditions and preparation of working solutions

Attention:

Buffers BQ1 and BW contain chaotropic salt! Wear gloves and goggles!

All components of the NucleoSpin® 96 DNA FFPE kit should be stored at room temperature (18–25 °C) and are stable for up to one year.

Before starting any NucleoSpin® DNA FFPE NucleoSpin® 96 DNA FFPE protocol, prepare the following:

Wash Buffer B5: Add the indicated volume of ethanol (96–100%) to **Buffer B5 Concentrate** before use. Mark the label of the bottle to indicate that ethanol was added. Store **Wash Buffer B5** at room temperature (18–25 °C) for up to one year.

Proteinase K: Before first use of the kit, add indicated volume of Proteinase Buffer PB to each vial of the **lyophilized Proteinase K**. Dissolved Proteinase K solution should be stored at -20 °C and is stable for 6 months

NucleoSpin® 96 DNA FFPE		
REF	1 x 96 preps 740240.1	4 x 96 preps 740240.4
Wash Buffer B5 (Concentrate)	100 mL Add 400 mL ethanol	2 x 100 mL Add 400 mL ethanol to each bottle
Proteinase K (lyophilized)	1 vial (75 mg) Add 2.8 mL Proteinase Buffer	4 vials (75 mg/vial) Add 2.8 mL Proteinase Buffer to each vial




4 Safety instructions

The following components of the **NucleoSpin® 96 DNA FFPE** kit contain hazardous chemicals.

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

GHS classification

Only harmful features need not be labeled with H and P phrases until 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
Inhalt	Gefahrstoff	GHS-Symbol	H-Sätze	P-Sätze
BQ1	Guanidine hydrochloride 50–66 % <i>Guanidinhydrochlorid 50–66 %</i> CAS 50-01-1	 WARNING ACHTUNG	302, 315, 319	280, 301+312, 302+352, 305+351+338, 330, 332+313, 337+313
BW	Guanidine hydrochloride 36–50 % + isopropanol 20–50 % Guanidinhydrochlorid 36–50 % + Isopropanol 20–50 % CAS 50-01-1, CAS 67-63-0	 WARNING ACHTUNG	226, 302, 319	210, 233, 280, 301+312, 305+351+338, 330, 337+313, 403+235
Proteinase K	Proteinase K, lyophilized <i>Proteinase K, lyophilisiert</i> CAS 39450-01-6	 DANGER GEFAHR	315, 319, 334, 335	261, 280, 302+352, 304+340, 305+351+338, 312, 332+313, 337+313, 342+311, 403+233

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 315	Causes skin irritation. <i>Verursacht Hautreizungen.</i>
H 319	Causes serious eye irritation. <i>Verursacht schwere Augenreizung.</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 335 May cause respiratory irritation.
Kann die Atemwege reizen.

Precaution phrases

- P 210 Keep away from heat, hot surfaces, sparks, open flames and other ignition sources. No smoking.
Von Hitze, heißen Oberflächen, Funken, offenen Flammen sowie anderen Zündquellenarten fernhalten. Nicht rauchen.
- P 233 Keep container tightly closed.
Behälter dicht verschlossen halten.
- P 261 Avoid breathing dust.
Einatmen von Staub vermeiden.
- P 280 Wear protective gloves / eye protection.
Schutzhandschuhe / Augenschutz tragen.
- P 301+312 IF SWALLOWED: Call a POISON CENTER / doctor / ... / if you feel unwell.
BEI VERSCHLUCKEN: Bei Unwohlsein GIFTINFORMATIONSZENTRUM / Arzt / ... anrufen.
- P 302+352 IF ON SKIN: Wash with plenty of water / ...
BEI KONTAKT MIT DER HAUT: Mit viel Wasser / ... waschen.
- P 304+340 IF INHALED: Remove victim to fresh air and keep at rest in a position comfortable for breathing.
BEI EINATMEN: An die frische Luft bringen und in einer Position ruhigstellen, die das Atmen erleichtert.
- P 305+351+338 IF IN EYES: Rinse continuously with water for several minutes. Remove contact lenses if present and easy to do – continue rinsing.
BEI KONTAKT MIT DEN AUGEN: Einige Minuten lang behutsam mit Wasser spülen. Vorhandene Kontaktlinsen nach Möglichkeit entfernen. Weiter spülen.
- P 312 Call a POISON CENTER / doctor / ... / if you feel unwell.
Bei Unwohlsein GIFTINFORMATIONSZENTRUM / Arzt / ... anrufen.
- P 330 Rinse mouth.
Mund ausspülen.
- P 332+313 If skin irritation occurs: Get medical advice / attention.
Bei Hautreizung: Ärztlichen Rat einholen / ärztliche Hilfe hinzuziehen.
- P 337+313 Get medical advice / attention.
Bei anhaltender Augenreizung: Ärztlichen Rat einholen / ärztliche Hilfe hinzuziehen.
- P 342+311 If experiencing respiratory symptoms: Call a POISON CENTER / doctor / ...
Bei Symptomen der Atemwege: GIFTINFORMATIONSZENTRUM / Arzt / ... anrufen.
- P 403+233 Store in a well ventilated place. Keep container tightly closed.
Behälter dicht geschlossen an einem gut belüfteten Ort aufbewahren.
- P 403+235 Store in a well ventilated place. Keep cool.
Kühl an einem gut belüfteten Ort aufbewahren.

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

6	Bind DNA to silica membrane of the NucleoSpin® FFPE DNA Binding Plate	5,600 x g, 10 min
7	Wash silica membrane	500 µL BW 5,600 x g, 2 min 700 µL B5 5,600 x g, 4 min
8	Dry silica membrane	70 °C, 10 min
9	Elute DNA	100 µL BE (70 °C) 5,600 x g, 2 min <i>Optional: Repeat elution step once.</i>

Detailed protocol

- For hardware requirements, refer to section 2.4.

Before starting the preparation:

- Check if Buffer B5 and Proteinase K were prepared according to section 3.
- Set incubator or oven to 60 °C.
- Preheat Elution Buffer BE to 70 °C.

1 Deparaffinize sample

Place each sample into a suitable 1.5 mL microreaction tube.

Add **400 µL Paraffin Dissolver** to the sample. Incubate **3 min** at **60 °C** (to melt the paraffin). **Vortex or shake** the sample immediately (at 60 °C) at a vigorous speed to dissolve the paraffin. Cool down sample to room temperature.

Make sure that paraffin completely melts during the heat incubation step and mix well after melting to completely dissolve the paraffin.

Insufficient mixing of the heated sample may cause recurrence of solid paraffin particles. Make sure the sample does not comprise more than 15 mg paraffin or adjust the volume of Paraffin Dissolver.

2 Lyse sample

Add **200 µL Lysis Buffer FL** and **25 µL Proteinase K** to the sample. Mix well by repeated pipetting up and down or pulse vortexing or shaking, centrifuge at **11,000 x g** for **1 min** and incubate at **56 °C** for **1–3 h** or overnight with shaking.

3 Decrosslink

Set heating block to 90 °C.

Centrifuge at **11000 x g** for **30 s**.

Decrosslinking step is strongly recommended for short time lysis (< 3 h) and may be omitted after overnight lysis.

Add **100 µL Buffer D-Link, mix** by repeated pipetting up and down, vortexing for 5 s or shaking.

Centrifuge for **1 min** at **11000 x g** to obtain phase formation.

Incubate at **90 °C** for **30 min**, **mix** by vortexing for 5 s. **Cool** down samples to room temperature. Centrifuge samples for **1 min** at **11000 x g**.

Transfer 300 µL of the lower aqueous phase from each sample to a Square-well Block for further processing.

4 Adjust DNA binding conditions

Add **300 µL Buffer BQ1** and **300 µL 96–100 % ethanol** to each sample. Take care not to moisten the rims of the individual wells while dispensing the buffer. Mix by pipetting up and down 5 times.

Ethanol and Buffer BQ1 can be premixed before addition to the samples, if the mixture is to be used during the next 3 months. Never centrifuge at higher g-forces or for longer periods as DNA will precipitate.

5 Transfer lysates

Place a NucleoSpin® DNA FFPE Binding Plate on a MN 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.

Transfer the lysates resulting from step 4 carefully into the wells of the NucleoSpin® FFPE DNA Binding Plate. Continue with the next samples. Do not moisten the rims of the individual wells while dispensing the samples – moistened rims may cause cross contamination during centrifugation. After transfer, seal the openings of the plate with Self-adhering PE Foil.

6 Bind DNA to silica membrane

Place the MN Square-well Blocks with NucleoSpin® FFPE DNA Binding Plates onto the centrifuge carriers and insert them into the rotor buckets. Centrifuge at **5,600–6,000 x g** for **2 min**.

Typically, the lysates will have passed through the silica membrane within a few minutes. The centrifugation process can be extended to 10 min, if the lysates have not passed completely.

7 Wash silica membrane

1st wash

Remove the Self-adhering PE Foil and add **500 µL Buffer BW** to each well of the NucleoSpin® FFPE DNA Binding Plate. Seal the plate with a new Self-adhering PE Foil and centrifuge again at **5,600–6,000 x g** for **2 min**.

2nd wash

Remove the Self-adhering PE Foil and add **700 µL Buffer B5** to each well of the NucleoSpin® FFPE DNA Plate. Seal the plate with a new Self-adhering PE Foil and centrifuge again at **5,600–6,000 x g** for **4 min**.

During this step, as much ethanolic Buffer B5 as possible is removed by centrifugation.

8 Dry silica membrane

Remove the Self-adhering PE Foil and place the NucleoSpin® FFPE DNA Binding Plate on an opened Rack of Tube Strips. Place it in an incubator for **10 min at 70 °C** to evaporate residual ethanol.

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

Note: The ethanol in Buffer B5 may inhibit enzymatic reactions and should be removed completely before eluting DNA.

9 Elute DNA

Dispense **100 µL** preheated **Buffer BE** (70 °C) to each well of the NucleoSpin® FFPE DNA Binding Plate. Dispense the buffer directly onto the membrane. Incubate at room temperature for 1 min.

Centrifuge at **5,600–6,000 x g** for **2 min**. *Optional: Repeat elution step once.*

Remove the NucleoSpin® FFPE DNA Binding Plate from the Rack of Tube Strips by lifting the plate at one side carefully as the Tube Strips may stick to the outlets of the NucleoSpin® FFPE DNA Binding Plate. For alternative elution procedures see section 2.6.

Alternatively, use the Round-well Block to collect the eluted DNA.

5.2 Vacuum processing

Protocol-at-a-glance

For hardware requirements refer to section 2.4.
For detailed information on each step see page 20.

Before starting the preparation:

- Check if Proteinase K and buffer B5 were prepared according to section 3.

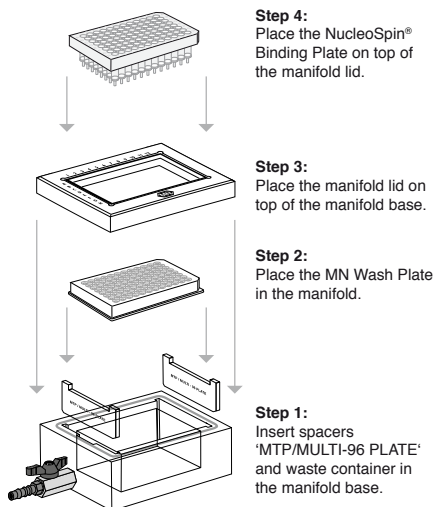
1	400 µL Paraffin Dissolver
Deparaffinize sample	60 °C, 3 min
	Mix
	Cool down to RT
2	Add to each sample
Lyse sample	200 µL FL
	25 µL Proteinase K
	Mix
	11,000 x g, 1 min
	56 °C, for 1–3 h or overnight
3	11,000 x g, 30 s
Decrosslink	100 µL D-Link
	mix
	11,000 x g, 30 s
	90 °C, 30 min
Load sample to processing plate	11,000 x g, 1 min
	Load aqueous phase (lower) in Square-well Block (approx. 300 µL)
4	300 µL BQ1
Adjust DNA binding conditions	300 µL Ethanol (96-100 %)
	Mix
	Prepare the NucleoVac 96 Vacuum Mantifold
5	Transfer lysates to NucleoSpin® DNA FFPE Binding plate

6	Bind DNA to silica membrane of the NucleoSpin® FFPE DNA Binding Plate	- 0.2 bar*, 5 min
7	Wash silica membrane	600 µL BW 900 µL B5 900 µL B5 - 0.2 bar*, 5 min each step Remove MN Wash Plate
8	Dry silica membrane	- 0.6 bar*, 10 min
9	Elute DNA	100 µL BE (70 °C) - 0.4 bar*, 2 min
<i>Optional: Repeat elution step once</i>		

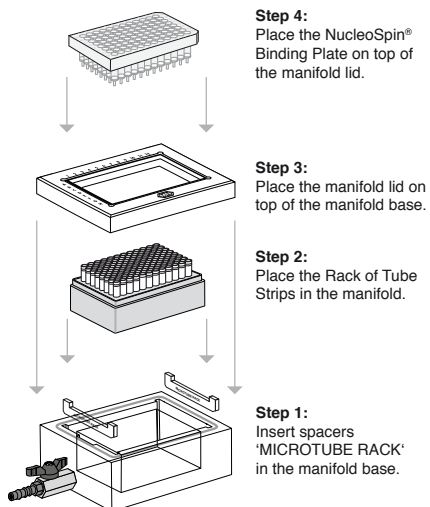
* Reduction of atmospheric pressure

Setup of vacuum manifold:

Binding / Washing steps



Elution step



Detailed protocol

- For hardware requirements, refer to section 2.4.

Before starting the preparation:

- Check if Buffer B5 and Proteinase K were prepared according to section 3.
- Set incubator or oven to 60 °C.
- Preheat Elution Buffer BE to 70 °C.

1 Deparaffinize sample

Place the sample into a suitable 1.5 mL microreaction tube or into a Round-well Block..

Add **400 µL Paraffin Dissolver** to the sample. Incubate **3 min at 60 °C** (to melt the paraffin). **Vortex or shake** the sample immediately (at 60 °C) at a vigorous speed to dissolve the paraffin. Cool down sample to room temperature.

Make sure that paraffin completely melts during the heat incubation step and mix well after melting to completely dissolve the paraffin. Insufficient mixing of the heated sample may cause recurrence of solid paraffin particles. Make sure the sample does not comprise more than 15 mg paraffin or adjust the volume of Paraffin Dissolver.

2 Lyse sample

Add **200 µL Lysis Buffer FL** and **25 µL Proteinase K** to the sample. Mix well by repeated pipetting up and down or pulse vortexing or shaking, centrifuge at **11,000 x g** for **1 min** and incubate at **56 °C** for **1–3 h** or overnight with shaking.

3 Decrosslink

Set heating block to 90 °C.

Centrifuge at **11000 x g** for **30 s**.

Decrosslinking step is strongly recommended for short time lysis (< 3 h) and may be omitted after overnight lysis.

Add **100 µL Buffer D-Link**, **mix** by repeated pipetting up and down, vortexing for 5 s or shaking.

Centrifuge for **1 min** at **11000 x g** to obtain phase formation.

Incubate at **90 °C** for **30 min**, **mix** by vortexing for 5 s. **Cool** down samples to room temperature. Centrifuge samples for **1 min** at **11000 x g**.

Transfer 300 µL of the lower aqueous phase from each sample to a Square-well Block for further processing.

4 Adjust DNA binding conditions

Add **300 µL Buffer BQ1** and **300 µL 96–100 % ethanol** to each sample. Take care not to moisten the rims of the individual wells while dispensing the buffer. Mix by pipetting up and down 5 times.

Ethanol and Buffer BQ1 can be premixed before addition to the samples, if the mixture is to be used during the next 3 months. Never centrifuge at higher g-forces or for longer periods as DNA will precipitate.

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® FFPE DNA Binding Plate on top of the manifold

5 Transfer lysates

Place a NucleoSpin® DNA FFPE Binding Plate on a MN 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.

Transfer the lysates resulting from step 4 carefully into the wells of the NucleoSpin® FFPE DNA Binding Plate. Continue with the next samples. Do not moisten the rims of the individual wells while dispensing the samples – moistened rims may cause cross contamination during centrifugation. After transfer, seal the openings of the plate with Self-adhering PE Foil.

6 Bind DNA to silica membrane

Apply vacuum until all lysates have passed through the wells of the NucleoSpin® FFPE DNA Binding Plate (**-0.2 bar***; **5 min**). Release the vacuum.

7 Wash silica membrane

1st wash

Add **600 µL Buffer BW** to each well of the NucleoSpin® FFPE DNA Binding Plate. Apply vacuum (**-0.2 bar***; **5 min**) until all buffer has passed through the wells of the NucleoSpin® FFPE DNA Binding Plate. Release the vacuum.

2nd wash

Add **900 µL Buffer B5** to each well of the NucleoSpin® FFPE DNA Binding Plate. Apply vacuum (**-0.2 bar***; **5 min**) until all buffer has passed through the wells of the NucleoSpin® FFPE DNA Binding Plate. Release the vacuum.

* Reduction of atmospheric pressure

3rd wash

Add **900 µL Buffer B5** to each well of the NucleoSpin® Tissue Binding Plate. Apply vacuum (**- 0.2 bar***; **5 min**) until all buffer has passed through the wells of the NucleoSpin® FFPE DNA Binding Plate. Release the vacuum.

Remove MN Wash Plate

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

8 Dry silica membrane

Insert the NucleoSpin® FFPE DNA Binding Plate into the lid, and close the manifold.

Apply maximum vacuum (**at least -0.6 bar***) for **10 min** to dry the membrane completely. This step is necessary to eliminate traces of ethanol.

Note: The ethanol in Buffer B5 inhibits enzymatic reactions and has to be removed completely before eluting DNA.

Finally, release the vacuum.

9 Elute DNA

Insert spacers “Microtube rack” into the NucleoVac Vacuum Manifold’s short sides. Place a Rack of Tube Strips onto the spacer. Close the vacuum manifold and place the NucleoSpin® FFPE DNA Binding Plate on top. Dispense **100 µL** preheated **Buffer BE** onto the membrane. Incubate for **3 min** at **room temperature**. Apply vacuum for elution (**- 0.4 bar***; **2 min**). Release the vacuum and repeat the elution step once.

For alternative elution procedures see section 2.6.

Finally, close the Tube Strips with Cap Strips for storage. Centrifuge Rack of Tube Strips briefly to collect all sample at the bottom of the Tube Strips.

Alternatively, use the Round-well Block for elution. Insert spacers “Round well Block” into the manifold.

6 Appendix

6.1 Troubleshooting

Problem	Possible cause and suggestions
No or poor DNA yield	<p><i>Incomplete lysis</i></p> <ul style="list-style-type: none">• 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 FL/Proteinase K mixture. Incubate until the samples are completely lysed. <p><i>Reagents not applied properly</i></p> <ul style="list-style-type: none">• Prepare Buffer B5 and Proteinase K solution according to instructions (see section 3). Add Buffer BQ1 and ethanol to the lysates before loading them to the wells of the NucleoSpin® FFPE DNA Binding Plate. <p><i>Suboptimal elution of DNA from the column</i></p> <ul style="list-style-type: none">• Preheat Buffer BE to 70 °C before elution. Apply Buffer BE 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 BE (pH 8.5).
RNA contamination	<p><i>RNA in sample</i></p> <ul style="list-style-type: none">• If RNA-free DNA 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, section 6.2). Incubate for 15 min with moderate shaking.

Poor performance of genomic DNA in enzymatic reactions	<p><i>Carry-over of ethanol</i></p> <ul style="list-style-type: none">• After washing with Buffer B5, centrifuge ≥ 4 min at 5,600–6,000 x <i>g</i> in order to remove ethanolic Buffer B5 completely and evaporate residual ethanol by incubating the NucleoSpin® DNA FFPE Binding Plate at 70 °C for 10 min.• Increase vacuum drying time to 15 min. <p><i>Contamination of DNA with inhibitory substances</i></p> <ul style="list-style-type: none">• Do not elute DNA with TE buffer. EDTA may inhibit enzymatic reactions. Repurify DNA and elute in Buffer BE.
Insufficient Paraffin removal	<p><i>Paraffin amount too high</i></p> <p>Using too many paraffin sections or excess of paraffin may cause immediate solidification of the paraffin after heat incubation when cooling lysate down to room temperature. Repeat heat incubation with increased (e.g., doubled) volume of paraffin dissolver.</p>

6.2 Ordering information

Product	REF	Pack of
NucleoSpin® 96 DNA FFPE	740240.1	1 x 96 preps
	740240.4	4 x 96 preps
NucleoMag® DNA FFPE	744320.1	1 x 96 preps
	744320.4	4 x 96 preps
NucleoSpin® FFPE DNA XS	740980.10/.50/.250	10/50/250 preps
Paraffin Dissolver (blue)	740343.60	60 mL
Decrosslink Buffer D-Link	740979.30	30 mL
Buffer BQ1	740923.1	1 L
Buffer B5 Concentrate (for 500 mL Buffer B5)	740921.100	100 mL
Buffer BW	740922.500	500 mL
Proteinase K (lyophilized)	740506	100 mg
RNase A (lyophilized)	740505.50	50 mg
	740505	100 mg
NucleoVac 96 Vacuum Manifold	740681	1
NucleoVac Vacuum Regulator	740641	1
Self-adhering PE Foil	740676	50

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

6.3 Product use restriction/warranty

NucleoSpin® 96 DNA FFPE 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

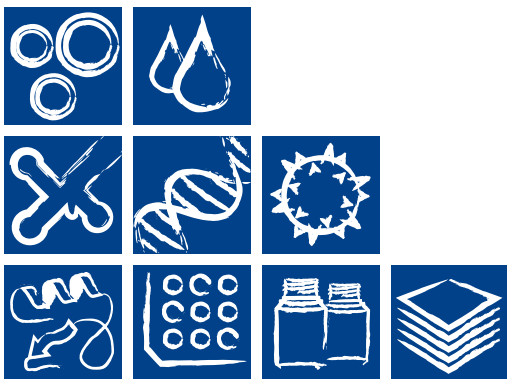
tech-bio@mn-net.com

Trademarks:

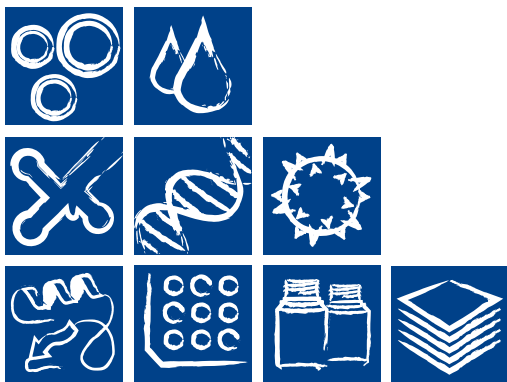
NucleoSpin is a registered trademark of MACHEREY-NAGEL GmbH & Co KG

PicoGreen is a registered trademark of Molecular Probs Inc.

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



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



MACHEREY-NAGEL



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

Germany

and international:

Tel.: +49 24 21 969-0

E-mail: info@mn-net.com

Switzerland:

MACHEREY-NAGEL AG

Tel.: +41 62 388 55 00

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

France:

MACHEREY-NAGEL EURL

Tel.: +33 388 68 22 68

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

USA:

MACHEREY-NAGEL Inc.

Tel.: +1 484 821 0984

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