



PCR clean-up

User manual

NucleoSpin[®] 96 PCR Clean-up

NucleoSpin[®] 96 PCR Clean-up Core Kit

This product distributed by Clontech Laboratories, Inc.

A Takara Bio Company

1290 Terra Bella Ave., Mountain View, CA 94043

www.clontech.com

For ordering information

1.800.662.2566

orders@clontech.com

For technical support

1.800.662.2566

tech@clontech.com



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

April 2014 / Rev. 04

MACHEREY-NAGEL

www.mn-net.com



Table of contents

1	Components	4
1.1	Kit contents	4
1.2	Reagents to be supplied by user	5
2	Product description	6
2.1	The basic principle	6
2.2	Kit specifications	6
2.3	Required hardware	8
2.4	Recommended accessories for use of the NucleoSpin® 96 PCR Clean-up Core Kit	8
2.5	Automated processing on robotic platforms	10
2.6	Elution procedures	10
3	Storage conditions and preparation of working solutions	12
4	Safety instructions	13
5	Protocols	14
5.1	NucleoSpin® 96 PCR Clean-up – manual vacuum processing	14
5.2	NucleoSpin® 96 PCR Clean-up – elution of DNA using a centrifuge	20
6	Appendix	21
6.1	Troubleshooting	21
6.2	Ordering information	22
6.3	References	23
6.4	Product use restriction/warranty	23

1 Components

1.1 Kit contents

NucleoSpin® 96 PCR Clean-up				
REF	1 x 96 preps	2 x 96 preps	4 x 96 preps	24 x 96 preps ¹
REF	740658.1	740658.2	740658.4	740658.24
Binding Buffer NT	25 mL	75 mL	2 x 75 mL	12 x 75 mL
Wash Buffer NT3 (Concentrate) ²	100 mL	2 x 100 mL	2 x 100 mL	12 x 100 mL
Elution Buffer NE ³	30 mL	60 mL	125 mL	6 x 125 mL
NucleoSpin® PCR Clean-up Binding Plate (yellow rings)	1	2	4	24
MN Wash Plate ⁴	1	2	4	24
Elution Plate U-bottom ⁵	1	2	4	24
User manual	1	1	1	6

¹ The kit for 24 x 96 preparations (REF 740658.24) consists of 6 x REF 740658.4.

² For preparation of working solutions and storage conditions see section 3.

³ Composition of Elution Buffer NE: 5 mM Tris/HCl, pH 8.5

⁴ Including six paper sheets

⁵ Including one Self-adhering PE foil

NucleoSpin® 96 PCR Clean-up Core Kit	
REF	4 x 96 preps 740464.4
Binding Buffer NT	2 x 75 mL
Wash Buffer NT3 (Concentrate) ¹	2 x 100 mL
Elution Buffer NE ²	125 mL
NucleoSpin® PCR Clean-up Binding Plate (yellow rings)	4
User manual	1

1.2 Reagents to be supplied by user

- 96–100 % ethanol

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

² Composition of Elution Buffer NE: 5 mM Tris/HCl, pH 8.5

2 Product description

2.1 The basic principle

The **NucleoSpin® 96 PCR Clean-up** kit allows direct clean-up of PCR reaction mixtures. Within the procedure the addition of chaotropic salt (Buffer NT) allows a reversible adsorption of the PCR products to the silica membrane of the NucleoSpin® 96 PCR Clean-up Binding Plates. High purity of the PCR-products is achieved by complete removal of primers, primer-dimers, salts, nucleotides, and proteins (e.g., polymerases, BSA) in subsequent washing steps using Buffer NT3. Highly pure PCR products are finally eluted with Elution Buffer NE (5 mM Tris/HCl, pH 8.5) or water (pH 8.5), and can be used directly for further applications.

2.2 Kit specifications

- **NucleoSpin® 96 PCR Clean-up** is designed for the fast 96-well purification of PCR products in the microtiter plate format (e.g., desalination, removal of enzymes, nucleotides and /or labeling reagents like biotin or radioactive ATP).
- The kit is for use with the NucleoVac 96 vacuum manifold (see ordering information) or similar suitable vacuum manifolds (see section 2.4).
- If using less than 96 samples the rubber pad or Self-adhering PE Foil (see ordering information) must be used in order to cover non used wells to maintain sufficient vacuum.
- The kit provides reagents and consumables for purification of up to 15 µg highly pure PCR products.
- Eluted PCR products are ready-to-use for several applications including automated fluorescent sequencing, cloning, or microarray technology.
- The **NucleoSpin® 96 PCR Clean-up** kit allows for the simultaneous processing of up to 96 samples typically within 45 minutes.

Kit specifications at a glance

Parameter	NucleoSpin® 96 PCR Clean-up
Format	96-well plates
Processing	Manual and automated, vacuum
Sample material	< 100 µL PCR reaction mixture
Fragment size	65 bp–10 kbp
Typical recovery	75–95 %
A_{260}/A_{280}	1.70–1.80
Elution volume	75–150 µL
Preparation time	45 min/plate
Binding capacity	15 µg

2.3 Required hardware

The **NucleoSpin® 96 PCR Clean-up** kits can be used **manually** with the NucleoVac 96 Vacuum Manifold (see ordering information).

This kit is intended for use under vacuum. A support protocol for elution under centrifugation is included (see section 5.2).

A support protocol for complete processing under centrifugation is available from our technical service (tech-bio@mn-net.com).

2.4 Recommended accessories for use of the NucleoSpin® 96 PCR Clean-up Core Kit

The **NucleoSpin® 96 PCR Clean-up Core Kit** provides buffers and NucleoSpin® Binding Plates only. Accessory plates (e.g., elution plates) are not provided with the core kit. The user can individually select additional consumables from a variety of suitable accessory plates according to his requirements for highest flexibility.

For use of **NucleoSpin® 96 PCR Clean-up Core Kit** follow the protocols (see section 5.1 or 5.2, respectively).

Recommended accessories for use of the **NucleoSpin® 96 PCR Clean-up Core Kit** are available from MACHEREY-NAGEL. For ordering information please refer to section 6.2.

2.5 Automated processing on robotic platforms

NucleoSpin® 96 PCR Clean-up can be used fully automated on many common laboratory workstations. For the availability of scripts and general considerations about adapting **NucleoSpin® 96 PCR Clean-up** on a certain workstation, please contact MN. Full processing under vacuum enables complete automation without the need for centrifugation steps, regarding drying of the binding membrane and elution step.

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® PCR Clean-up Binding Plate.

Drying of the NucleoSpin® PCR Clean-up Binding Plate under vacuum is sufficient because the bottom of the plate is protected from residues of wash buffer during the washing steps by the MN Wash Plate. As a result, we recommend trying to integrate the MN Wash Plate into the automated procedure. The MN Frame (see ordering information) 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 DNA. 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. Several application notes of the **NucleoSpin® 96 PCR Clean-up** kit (formerly known as NucleoSpin® 96 Extract II) on various liquid handling instruments can also be found at www.mn-net.com at Bioanalysis/Literature.

2.6 Elution procedures

Elution of purified PCR products: The efficiency of the DNA elution depends on the pH of the elution buffer. Elution is most effective at pH 8.0–8.5. When using nuclease-free water for elution, the pH value should be checked and, if necessary, adjusted to 8.0–8.5. Lower pH of the selected elution buffer may lead to lower recoveries. Yield of larger DNA fragments (> 5–10 kbp) can be increased by using pre-warmed (70 °C) elution buffer (also see Table below). An elution volume of 75–125 µL Buffer NE, as well as a 3–5 min incubation at room temperature (18–25 °C) of the elution buffer on the silica membrane are recommended.

See the following table for correlation between the dispensed elution buffer volume and typical recoveries following the standard protocol.

The recommended dispense volume of elution buffer is 125 µL.

Correlation between dispensed elution buffer volume and typical recovery

Dispensed elution buffer	75 µL	100 µL	125 µL	150 µL	175 µL
Recovered elution buffer containing PCR-products	30 ± 5 µL	55 ± 5 µL	80 ± 5 µL	105 ± 5 µL	130 ± 5 µL

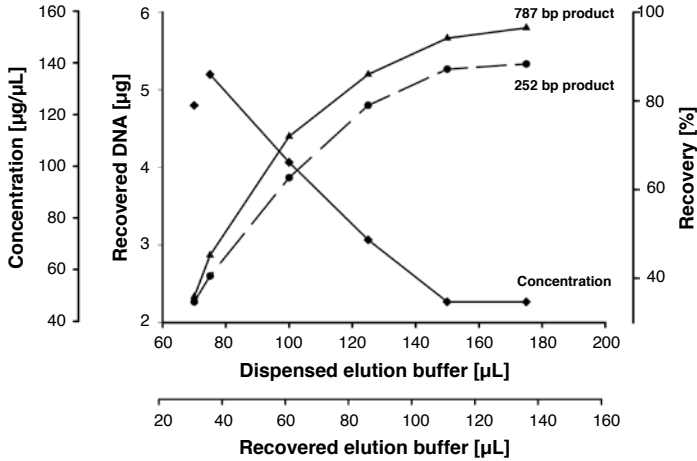


Figure 1: Recovery rate and concentration depend on elution volume.

Two different PCR products (252 bp and 787 bp) have been purified with the NucleoSpin® 96 PCR Clean-up kit.

Average DNA recovery rate depends on the size of PCR product	
Size of PCR product	Average DNA recovery rate
64 bp	60–80 %
164 bp	70–85 %
200 bp	70–85 %
490 bp	85–95 %
982 bp	85–95 %
1500 bp	80 %
2000 bp	75 %
4000 bp	50–60 %

3 Storage conditions and preparation of working solutions

Attention:

Storage conditions:

- **NucleoSpin® 96 PCR Clean-up/96 PCR Clean-up Core Kits** should be stored at room temperature (18–25 °C) and are stable for up to one year.

Before starting any **NucleoSpin® 96 PCR Clean-up/96 PCR Clean-up Core Kit** purification prepare the following:

- **Wash Buffer NT3:** Add the indicated volume of ethanol (96–100%) to **Buffer NT3 Concentrate**. Mark the label of the bottle to indicate that ethanol was added. Wash Buffer NT3 is stable at room temperature (18–25 °C) for up to one year.

NucleoSpin® 96 PCR Clean-up				
REF	1 x 96 preps 740658.1	4 x 96 preps 740658.2	4 x 96 preps 740658.4	24 x 96 preps 740658.24
Wash Buffer NT3 (Concentrate)	100 mL	2 x 100 mL	2 x 100 mL	12 x 100 mL
	Add 400 mL ethanol	Add 400 mL ethanol to each bottle	Add 400 mL ethanol to each bottle	Add 400 mL ethanol to each bottle

NucleoSpin® 96 PCR Clean-up Core Kit	
REF	4 x 96 preps 740464.4
Wash Buffer NT3 (Concentrate)	2 x 100 mL Add 400 mL ethanol


4 Safety instructions

The following components of the **NucleoSpin® 96 PCR Clean-up** and the **NucleoSpin® 96 PCR Clean-up Core Kits** contains hazardous contents. *Wear gloves and goggles and follow the safety instructions given in this section.*

GHS classification

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

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

Component	Hazard contents	GHS symbol	Hazard phrases	Precaution phrases
<i>Inhalt</i>	<i>Gefahrstoff</i>	<i>GHS Symbol</i>	<i>H-Sätze</i>	<i>P-Sätze</i>
NT	Guanidinium thiocyanate 30–60 % <i>Guanidiniumthiocyanat 30–60 %</i>	 Warning <i>Achtung</i>	302, 412, 031	260, 273, 301+312, 330

Hazard phrases

- H 031 Contact with acids liberates toxic gas.
Entwickelt bei Berührung mit Säure giftige Gase.
- H 302 Harmful if swallowed.
Giftig bei Verschlucken.
- H 412 Harmful to aquatic life with long lasting effects.
Schädlich für Wasserorganismen, mit langfristiger Wirkung.

Precaution phrases

- P 260 Do not breathe dust / fume / gas / mist / vapours / spray.
Staub / Rauch / Gas / Nebel / Dampf / Aerosol nicht einatmen.
- P 273 Avoid release to the environment.
Freisetzung in die Umwelt vermeiden. Mund ausspülen.
- P 301+312 IF SWALLOWED: Call a POISON CENTER/ doctor/.../if you feel unwell.
Bei Verschlucken: Bei Unwohlsein GIFTINFORMATIONSZENTRUM / Arzt /... anrufen.
- P 330 Rinse mouth.
Mund ausspülen.

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

5 Protocols

5.1 NucleoSpin® 96 PCR Clean-up – manual vacuum processing

- For hardware requirements refer to section 2.3.
- For detailed information regarding the vacuum manifold set-up see page 17.
- For detailed information on each step see page 18.
- For use of the NucleoSpin® 96 PCR Clean-up [Core Kit](#) (REF 740464.4), refer to section 2.4 regarding recommended accessories.

Before starting the preparation:

- Check if Buffer NT3 was prepared according to section 3.
- Set up the vacuum manifold according to the scheme

Protocol-at-a-glance

1	Adjust the volume of the reaction mixture to 100 µL using Tris buffer (pH 7.0–7.5), nuclease-free water (pH 7.0–7.5), or use Buffer NE	For PCR samples < 100 µL
2	Dispense binding buffer to NucleoSpin® PCR Clean-up Binding Plate	200 µL NT
3	Transfer PCR samples to NucleoSpin® PCR Clean-up Binding Plate and mix	100 µL diluted PCR sample
4	Bind DNA to silica membrane of the NucleoSpin® PCR Clean-up Binding Plate by applying vacuum	-0.2 bar to -0.4 bar* (1 min)
5	Wash silica membrane	2 x 900 µL NT3 -0.2 bar to -0.4 bar* (1 min)
6	Remove MN Wash Plate	

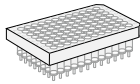
* Reduction of atmospheric pressure

7	Dry NucleoSpin® PCR Clean-up Binding Plate by applying vacuum <i>Optional: Dry the outlets of the NucleoSpin® PCR Clean-up Binding Plate by placing it on a sheet of filter paper before applying vacuum</i>	-0.3 to -0.4 bar* 10–15 min (run pump continuously)*
8	Insert Elution Plate U-bottom	
9	Elute DNA <i>Optional: Incubate 1–3 min</i>	75–150 µL NE -0.4 to -0.6 bar* (1 min)

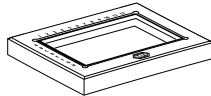
* Reduction of atmospheric pressure

Setup of vacuum manifold:

Binding / Washing / Elution steps



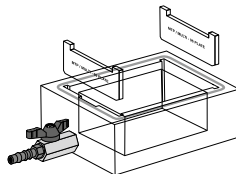
NucleoSpin® Binding Plate



MN Wash Plate



Elution Plate



Manifold base with spacers
'MTP/Multi-96 Plate' inserted

Binding / Washing step

Elution step

* Reduction of atmospheric pressure

Detailed protocol

For processing of NucleoSpin® 96 PCR Clean-up under vacuum the NucleoVac 96 Vacuum Manifold is required.

Before starting the preparation:

- Check if Buffer NT3 was prepared according to section 3.
-

1 Adjust the volume of reaction mixture

For PCR reaction volumes below 100 µL: Before starting the purification procedure, add Tris buffer (10 mM, pH 7.0), **nuclease-free water** (pH 7.0–7.5), or **Elution Buffer NE** to adjust the reaction mixture to a final volume of 100 µL.

Note: If less than 100 µL of PCR reaction mixture is used the volume of Binding Buffer NT has to be adjusted. It is mandatory that the ratio of Buffer NT : PCR reaction mixture is 2 : 1.

Prepare the NucleoVac 96 Vacuum Manifold

Insert spacers labeled 'MTP/Multi-96 Plate' notched side up into the grooves located on the short sides of the manifold. Insert waste container into manifold base. Place the MN Wash Plate on top of the spacers. Insert NucleoSpin® PCR Clean-up Binding Plate into the manifold lid and place lid on the manifold base. Close manifold base with the manifold lid. Close the vacuum manifold's valve, check and adjust the vacuum (-0.2 bar*).

2 Dispense binding buffer to the NucleoSpin® PCR Clean-up Binding Plate (column wise processing is recommended)

Add **200 µL Buffer NT** to each well of the NucleoSpin® PCR Clean-up Binding Plate.

3 Transfer PCR samples to the NucleoSpin® PCR Clean-up Binding Plate and mix

Mix by pipetting up and down 5 times.

Optional: Pre-mix PCR reaction and Buffer NT in a Square-well Block etc. (not supplied).

4 Bind DNA to silica membrane

Apply vacuum by opening the valve and press down the plate slightly until flow-through starts. Allow the samples to pass the columns and release vacuum by closing the valve **(-0.2 to -0.4 bar*, 1 min)**.

* Reduction of atmospheric pressure

5 Wash silica membrane

1st wash

Add 900 µL Buffer NT3 (with ethanol added) to each well of the NucleoSpin® PCR Clean-up Binding Plate.

Apply vacuum by opening the valve. Press down the NucleoSpin® PCR Clean-up Binding Plate slightly until flow-through starts. Allow the buffer to pass the columns. Release the vacuum (**-0.2 to -0.4 bar***, **1 min**).

2nd wash

Repeat this washing step once.

6 Remove MN Wash Plate

After the final washing step, close the valve, release the vacuum, and remove the NucleoSpin® PCR Clean-up Binding Plate. Remove manifold lid, MN Wash Plate, and waste container from the vacuum manifold.

7 Dry NucleoSpin® PCR Clean-up Binding Plate

Remove any residual washing buffer from the NucleoSpin® PCR Clean-up Binding Plate. If necessary, tap the outlets of the NucleoSpin® PCR Clean-up Binding Plate onto a clean paper sheet (supplied with the MN Wash Plate) or soft tissue until no drops come out. Insert the NucleoSpin® PCR Clean-up Binding Plate into the lid and close the manifold. Apply vacuum of **-0.3 to -0.4 bar*** for **at least 10 min** to dry the membrane completely. This step is necessary to eliminate traces of ethanol.

Note: The ethanol in Buffer NT3 inhibits enzymatic reactions and has to be completely removed before eluting the DNA.

Finally, close the release the vacuum.

* Reduction of atmospheric pressure

8 Insert Elution Plate U-bottom

Insert the Elution Plate U-bottom on the spacers inside the manifold base. For elution into microtiter plates spacers '*MTP/Multi-96 Plate*' are required which are already inserted into the manifold base from the previous steps. Reassemble the vacuum manifold as described before.

Or

Elution into Rack of Tube Strips (not provided with the kit, see ordering information): Insert spacers '*Microtube rack*', notched side up, into the grooves located at the short sides of the vacuum manifold. Rest the Rack of Tube Strips on the spacers inside the manifold base and reassemble the vacuum manifold as described before.

9 Elute DNA

Add **75–150 µL Elution Buffer NE** (5 mM Tris-HCl, pH 8.5) **or water** (pH 8.5) to each well of the NucleoSpin® PCR Clean-up Binding Plate.

The buffer should be dispensed onto the center of the silica membrane. Incubate for **1–3 min at room temperature** (optionally), apply vacuum, and collect the eluted DNA. After the elution buffer has passed the columns, release the vacuum (**-0.4 to -0.6 bar***, **1–2 min**).

Remove the Elution Plate U-bottom (or Rack of Tube Strips) containing eluted DNA and seal them with the supplied adhesive cover foil (or Cap Strips for Tube Strips) for further storage.

* Reduction of atmospheric pressure

5.2 NucleoSpin® 96 PCR Clean-up – elution of DNA using a centrifuge

Eluting the purified DNA in a centrifuge may be necessary when higher concentrations of the final DNA are required for downstream applications. Using a centrifuge allows the dispensed volume to be reduced down to 50–75 µL, giving a DNA concentration of about 70–200 ng/µL (depending on elution buffer volume and fragment length).

Required hardware:

- For centrifugation, a microtiterplate centrifuge that can accommodate the NucleoSpin® PCR Clean-up Binding Plate stacked on a rack of Tube Strips is required. It is also necessary that the centrifuge reached accelerations of 5,600–6,000 x *g* (bucket height: 85 mm).
 - For centrifugation, a microtiterplate centrifuge is required which is able to accommodate the NucleoSpin® PCR Clean-up Binding Plate stacked on Rack of Tube Strips and reaches accelerations of 5,600–6,000 x *g* (bucket height: 85 mm).
 - Suitable elution tubes: Rack of Tube Strips has to be ordered separately (see ordering information).
-

1 Stop the method after the final washing step with Buffer NT3. Remove the NucleoSpin® PCR Clean-up Binding Plate from the manifold's top and tap on a sheet of filter paper to remove residual wash buffer from the outlets.

2 Place the NucleoSpin® PCR Clean-up Binding Plate on top of a MN Square-well Block (not supplied with the kit, see ordering information) and centrifuge for **10 min** at **maximum speed** (> 4,000 x *g*, optimal 5,800 x *g*).

*Note: Do not use a microtiter plate as a support for the NucleoSpin® PCR Clean-up Binding Plate. Microtiter plates may crack under centrifugation at > 1,500 x *g*.*

3 Place the NucleoSpin® PCR Clean-up Binding Plate on top of a Rack of Tube Strips (not supplied with the kit, see ordering information). Dispense **Elution Buffer NE** (50–150 µL) directly onto the silica membrane and incubate for **1–3 min** at **room temperature**.

4 Centrifuge for **2 min** at **maximum speed** (> 4,000 x *g*, optimal 5,800 x *g*) to collect the DNA.

Remove the Rack of Tube Strips containing eluted DNA and close them with Cap Strips for further storage.

6 Appendix

6.1 Troubleshooting

Problem	Possible cause and suggestions
Poor DNA yield	<p data-bbox="348 344 981 392"><i>No ethanol added to Buffer NT3 Concentrate, ethanol evaporated</i></p> <ul data-bbox="348 405 981 481" style="list-style-type: none"> <li data-bbox="348 405 981 481">• Add indicated volume of ethanol to Buffer NT3 Concentrate and mix. Keep bottle tightly closed to prevent evaporation of ethanol. <p data-bbox="348 523 684 545"><i>Elution conditions are not optimal</i></p> <ul data-bbox="348 558 981 663" style="list-style-type: none"> <li data-bbox="348 558 981 663">• If possible, use a slightly alkaline elution buffer like Buffer NE (5 mM Tris-HCl, pH 8.5). When using nuclease-free water for elution, make sure the pH value is 8.5. Elution efficiencies drop dramatically at pH < 7. <p data-bbox="348 705 698 727"><i>Elution buffer volume is insufficient</i></p> <ul data-bbox="348 740 981 788" style="list-style-type: none"> <li data-bbox="348 740 981 788">• Optimal elution is achieved for an elution buffer volume of 100–150 μL. Do not use less than 75 μL elution buffer.
Suboptimal performance of PCR product in sequencing reactions, problems with downstream applications	<p data-bbox="348 826 564 849"><i>Carry-over of ethanol</i></p> <ul data-bbox="348 861 981 938" style="list-style-type: none"> <li data-bbox="348 861 981 938">• Be sure to remove all of ethanolic Buffer NT3 after the final washing step. Dry the NucleoSpin[®] PCR Clean-up Binding Plate for at least 10 min with maximum vacuum. <p data-bbox="348 979 736 1002"><i>Elution of PCR products with TE buffer</i></p> <ul data-bbox="348 1015 981 1145" style="list-style-type: none"> <li data-bbox="348 1015 981 1145">• EDTA may inhibit enzymatic reactions like DNA sequencing. Repurify the PCR products and elute with NE buffer or nuclease-free water. Alternatively, the DNA may be precipitated with ethanol and redissolved in buffer NE buffer or nuclease-free water. <p data-bbox="348 1187 818 1209"><i>Not enough DNA used in sequencing reactions</i></p> <ul data-bbox="348 1222 981 1273" style="list-style-type: none"> <li data-bbox="348 1222 981 1273">• Quantitate DNA by agarose gel electrophoresis before setting up sequencing reactions.

Problem	Possible cause and suggestions
Suboptimal performance of PCR product in sequencing reactions, problems with downstream applications (continued)	<i>Contamination of PCR product preparation with ethanol</i> <ul style="list-style-type: none"> Insufficient drying after final washing step with Buffer NT3. Remaining ethanol may cause problems with downstream applications like DNA sequencing or loading of samples onto agarose gel.
	<i>Eluted DNA contains residual primers/primer dimers</i> <ul style="list-style-type: none"> Minimize amount of primers in PCR reaction mixture. Make sure that the ratio of binding buffer NT:PCR reaction mixture is 2:1.

6.2 Ordering information

Product	REF	Pack of
NucleoSpin® 96 PCR Clean-up	740658 .1	1 x 96 preps
	740658 .2	2 x 96 preps
	740658 .4	4 x 96 preps
	740658 .24	24 x 96preps
NucleoSpin® 96 PCR Clean-up Core Kit	740464 .4	4 x 96 preps
NucleoSpin® 8 PCR Clean-up	740668	12 x 8 preps
	740668 .5	60 x 8 preps
NucleoSpin® 8 PCR Clean-up Core Kit	740463 .4	48 x 8 preps
MN Wash Plate	740479	4
	740479.24	24
Round-well Block with Cap Strips (set consists of 1 Round-well Block and 12 Cap Strips)	740475	4 sets
	740475.24	24 sets
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
Cap Strips	740478	48
	740478 .24	288
MN Square-well Block	740476	4
	740476 .24	24
Round-well Block Low (set consists of Round-well Block Low and Self-adhering Foil)	740487	4 sets
	740487.24	24 sets

Product	REF	Pack of
Elution Plate U-bottom (with Self-adhering Foil)	740486 .24	24 sets
Cap Strips	740638	30
Self-adhering PE Foil	740676	50
MN Frame	740680	1
NucleoVac 96 Vacuum Manifold	740681	1
NucleoVac Vacuum Regulator	740641	1

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

6.3 References

Vogelstein, B. & Gillespie, D. (1979) Proc. Natl. Acad. Sci. USA 76, 615-619.

6.4 Product use restriction/warranty

NucleoSpin® 96 PCR Clean-up (Core 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