

MACHEREY-NAGEL

User manual

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. ©2021



PCR clean-up Gel extraction

■ NucleoSpin® Gel and PCR Clean-up XS

February 2021 / Rev. 01

Table of contents

1 Components	4
1.1 Kit contents	4
1.2 Reagents, consumables, and equipment 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 Elution procedure	5
3 Storage conditions and preparation of working solutions	6
4 Safety instructions	7
5 Protocols	8
6 Appendix	10

1 Components

1.1 Kit contents

NucleoSpin® Gel and PCR Clean-up XS			
REF	10 preps 740611.10	50 preps 740611.50	250 preps 740611.250
Binding Buffer NTI	10 mL	40 mL	200 mL
Wash Buffer NT3 (Concentrate)*	6 mL	12 mL	50 mL
Elution Buffer NE**	13 mL	13 mL	13 mL
NucleoSpin® Gel and PCR Clean-up XS columns (yellow rings)	10	50	250
Collection Tubes (2 mL)	10	50	250
Short protocol	1	1	1

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

Reagents

- 96–100 % ethanol

Consumables

- 1.5 mL microcentrifuge tubes
- Disposable pipette tips

Equipment

- Manual pipettors
- Centrifuge for microcentrifuge tubes
- Heating block, water bath, or thermomixer for gel extraction
- Scalpel to cut agarose gels
- Vortex mixer
- Personal protection equipment (lab coat, gloves, goggles)

1.3 About this user manual

It is strongly recommended reading the detailed protocol sections of this user manual if the **NucleoSpin® Gel and PCR Clean-up XS** kit is used for the first time. Experienced users, however, may refer to the Protocol at a glance instead. All technical literature is available on the internet at www.mn-net.com.

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

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

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

2 Product description

2.1 The basic principle

NucleoSpin® Gel and PCR Clean-up XS is designed as a 2-in-1 kit to isolate DNA from enzymatic reactions or agarose gels. Binding buffer NT1 inactivates enzymes and promotes the binding of DNA to the silica layers inside the binding columns. Furthermore, Buffer NT1 is used in combination with an incubation at 50°C to melt agarose gels during DNA gel extraction. Bound DNA is stringently washed with buffer NT3 and eluted in elution volumes as low as 6 µl of Buffer NE, resulting in highly pure and concentrated DNA.

2.2 Kit specifications

- The NucleoSpin® Gel and PCR Clean-up XS kit is designed for research use only.

Table 1: Kit specifications at a glance

Parameter	NucleoSpin® Gel and PCR Clean-up XS
Format	Minispin, XS format
Processing	Manual
Sample material	≤ 200 µL enzymatic reaction ≤ 200 mg agarose gel
Elution volume	≥ 6 µL
Preparation time	< 15 min / 12 preps (enzymatic reaction clean up) 30 min / 12 preps (gel extraction)
Binding capacity	5 µg

2.3 Elution procedure

The XS column format is designed for low elution volumes, resulting in high DNA concentrations. Nevertheless there are tips and tricks to enhance the DNA recovery, especially when working with elution volumes below 10 µL.

The most relevant improvements in terms of nucleic acid recovery can be achieved by a two-fold elution. For a two-fold elution, the eluate of the elution step is reloaded onto the silica for a second elution procedure. Doing so, the total elution volume remains low while the recovery is enhanced.

This is especially true for gel extraction procedures. DNA tends to get stuck to the silica matrix and shows a delayed elution profile. Multiple rounds of elution can increase the DNA recovery significantly.

Keep in mind that chaotropic salt carry-over will increase the absorption at 230 nm without negative effects on downstream applications. Even minute amounts of thiocyanate, well below an influential concentration, show a strong absorption at 230 nm. There is no effect on the absorption at 260 nm.

3 Storage conditions and preparation of working solutions

Attention: Buffer NT1 contains guanidine thiocyanate! Wear gloves and goggles!

CAUTION: Buffer NT1 contains guanidine thiocyanate which can form highly reactive compounds when combined with bleach (sodium hypochloride). DO NOT add bleach or acidic solutions directly to the sample preparation waste.

- All kit components can be stored at room temperature (18–25°C) and are stable for at least one year.

Before starting the NucleoSpin® Gel and PCR Clean-up XS protocol prepare the following:


NucleoSpin® Gel and PCR Clean-up XS			
REF	10 preps 740611.10	50 preps 740611.50	250 preps 740611.250
Wash Buffer NT3 (Concentrate)	6 mL Add 24 mL ethanol	12 mL Add 48 mL ethanol	50 mL Add 200 mL ethanol

4 Safety instructions

The following components of the **NucleoSpin® Gel and PCR Clean-up XS** kits contain hazardous contents.

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

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
Inhalt	Gefahrstoff	GHS-Symbol	H-Sätze	P-Sätze
NTI	Guanidinium thiocyanate 30–45 % Guanidinthiocyanat 30–45 % CAS 593-84-0	 WARNING ACHTUNG	302, 412	264W, 273, 301+312, 330

Hazard phrases

- H 302 Harmful if swallowed.
Gesundheitsschädlich bei Verschlucken.
- H 412 Harmful in contact with skin.
Schädlich für Wasserorganismen, mit langfristiger Wirkung.

Precaution phrases

- P 264W Wash with water thoroughly after handling.
Nach Gebrauch mit Wasser gründlich waschen.
- P 273 Avoid release to the environment.
Freisetzung in die Umwelt vermeiden.
- P 301+312 IF SWALLOWED: Call a POISON CENTER / doctor / ... / if you feel unwell.
BEI VERSCHLUCKEN: Bei Unwohlsein GIFTINFORMATIONSZENTRUM/Arzt/... anrufen. Besondere Behandlung (siehe ... auf diesem Kennzeichnungsetikett).
- 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).



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

Before starting the preparation:

- Check if Wash Buffer NT3 was prepared according to section 3.
- When performing a gel extraction, excise DNA band from agarose gel and determine the weight of the gel slice. Preheat a heating block, thermomixer or water bath to 50°C.

1 Adjust DNA binding condition

Mix **1 volume of sample** with **2 volumes of Buffer NT1** (e. g., mix 20 µL PCR reaction with 40 µL NT1 or 150 mg agarose gel slice with 300 µL NT1).



**+ 2 vol NT1 per
1 vol sample**

Vortex

Agarose gel extraction only:

Incubate mixture at 50°C with constant shaking or repeated vortexing every 2–3 min until the gel is completely dissolved.

2 Bind DNA

Place a **NucleoSpin® Gel and PCR Clean-up XS Column** into a Collection Tube (2 mL) and load up to 500 µL sample.



Load sample

Centrifuge for **30 s** at **11,000 x g**. Discard flowthrough if it exceeds 100 µL and place the column back into the collection tube.



**11,000 x g
30 s**

If necessary, load remaining sample and repeat the centrifugation step.

3 Wash silica

Add **500 µL Buffer NT3** to the NucleoSpin® Gel and PCR Clean-up XS Column. Centrifuge for **30 s** at **11,000 x g**. Discard flowthrough and place the column into a 1.5 mL microcentrifuge tube (not supplied)



+ 500 µL NT3



**11,000 x g
30 sec**

4 Wash and dry silica

Add **300 µL Buffer NT3** to the NucleoSpin® Gel and PCR Clean-up XS column.



+ 300 µL NT3

Centrifuge for **1 min** at **11,000 x g**. Discard flowthrough and place the column into a 1.5 mL microcentrifuge tube (not supplied).



**11,000 x g
1 min**

5 Elute DNA

Add **6–12 µL Buffer NE** and incubate at room temperature (18–25°C) for **1 min**.

Centrifuge for **1 min at 11,000 x g**.

Reload the eluates onto the column and repeat the elution step.



+ 6–12 µL NE

**RT
1 min**

**11,000 x g
1 min**

6 Appendix

6.1 Troubleshooting

Problem	Possible cause and suggestions
Incomplete dissolving of gel slice	<p><i>Time and temperature</i></p> <ul style="list-style-type: none"> • Increase incubation time and vortex mixture repeatedly every 2–3 min until the gel slice is completely dissolved. Do not increase the incubation temperature to prevent loss of thermally denatured single stranded DNA
Low DNA yield	<p><i>Reagents prepared or stored improperly</i></p> <ul style="list-style-type: none"> • Add the indicated volume of 96–100 % ethanol to Buffer NT3. Mix well and keep bottle closed tightly. Prevent conditions where ethanol might evaporate (e. g., high temperatures, open bottles). <p><i>Incompletely dissolved gel slice</i></p> <ul style="list-style-type: none"> • Weigh agarose block and add 2 volumes of Buffer NT1. Vortex mixture continuously or every 2–3 min while melting the agarose gel. If necessary, increase incubation times, but do not increase incubation temperature. <p><i>Ethanol carry-over</i></p> <ul style="list-style-type: none"> • Take care that the columns' outlets or walls do not come in contact with the NT3 flowthrough after the second washing step. If in doubt, place the column back into the empty collection tube and re-centrifuge.

Problem	Possible cause and suggestions
Low DNA yield	<p><i>Incomplete elution</i></p> <ul style="list-style-type: none"> If DNA recovery is low, the remaining DNA is presumably still sticking to the silica. Losses during the binding or washing step are unlikely if the buffers are used, stored and prepared according to the protocol. To enhance the DNA recovery, it is recommended to either increase the total elution volume or to use repeated elution steps with the same aliquot, especially when using low elution volumes below 10 µL. Simply reload the eluate onto the silica and repeat the centrifugation.
	<p><i>Input quantification overestimated</i></p> <ul style="list-style-type: none"> Keep in mind, that photometry will sum up the absorption of any molecule which absorbs at a certain wavelength, e. g., 260 nm. A PCR reaction will contain not only the desired DNA amplicates, but also residual primers, unused dNTPs, enzymes, certain detergents or buffer substances, which all absorb at 260 nm and increase the measured "DNA" concentration. A photometric determination of the product recovery is not possible. To assess the product recovery, alternative techniques like gel electrophoresis, capillary electrophoresis or dsDNA specific fluorometry must be used. Depending on the PCR mix, 80–90 % of the initial absorption at 260 nm will be removed by the purification process.

Problem	Possible cause and suggestions
Low DNA yield	<p><i>Smear bands</i></p> <ul style="list-style-type: none"> Agarose gels might show artifacts when the agarose mesh is irregular due to insufficient homogenization of agarose in TAE /TBE buffer, differences in solidification, local thickness or other influences which alter the migration speed. The DNA itself might also be able to migrate in different conformations, depending on sequence and buffer temperature. While the main band might show a strong signal it is still possible that some fragments will migrate slightly faster or slower in an insufficient amount to give a detectable signal or show a smear. Excising the band closely will remove those fragments but is recommended nevertheless to reduce the amount of agarose.
Low $A_{260/230}$	<p><i>Chaotropic salt carry-over</i></p> <ul style="list-style-type: none"> To melt agarose gels and to bind nucleic acids to a silica surface, thiocyanate salts have proven effective and reliable. They have a long term track record showing successful purification of nucleic acids from various sample types and for various downstream applications. Nevertheless, these chaotropic salts express a very high absorption at 230 nm even at negligible concentrations as low as the micro molar range, while the absorption at 260 nm is zero. Thiocyanate salts will therefore bias the A_{260}/A_{230} purity ratio but not the DNA quantification. A low A_{260}/A_{230} is no indicator for suboptimal performance in downstream applications in this case and can be ignored. To improve the A_{260}/A_{230}, further washing steps with buffer NT3 can be performed. Additional Buffer NT3 will be necessary and can be ordered at MACHEREY-NAGEL (cat. no. 740598).

Problem	Possible cause and suggestions
High A_{320}	<p data-bbox="568 209 708 229"><i>Silica abrasion</i></p> <ul style="list-style-type: none"> <li data-bbox="568 244 969 469">• In rare cases, sheared silica fibers might pass the filter frit of the columns and are eluted. When measured in a photometer, those fibers will scatter the light, resulting in an overall overestimation of the absorption values. A high A_{320} value above 0.05 will be an indicator for particles which scatter light. <li data-bbox="568 483 975 710">• Silica abrasion can be easily pelleted by centrifugation and is usually already spun down by the elution. Vortexing and mixing the eluates might resuspend these particles again. To prevent this it is recommended to transfer the clear eluate without the easily visible white pellet into a new centrifuge tube.

Problem	Possible cause and suggestions
Suboptimal performance in downstream applications	<p><i>DNA was damaged by UV light</i></p> <ul style="list-style-type: none"> • Reduce UV exposure time to a minimum when excising a DNA fragment from an agarose gel <p><i>Ethanol carry-over</i></p> <ul style="list-style-type: none"> • Due to the low elution volumes, ethanol carry-over from the second washing step can reach a high concentration in the eluates and might inactivate enzymes. Take care that the outlets or walls of the columns do not come in contact with the NT3 flowthrough from the second washing step. If in doubt, discard flowthrough and place the columns back into the empty collection tubes. Repeat the centrifugation step with empty columns. <p><i>Silica abrasion</i></p> <ul style="list-style-type: none"> • Silica debris might scatter or quench the signals during photometric or fluorescence based measurements. Furthermore, the channels of capillary electrophoresis might be blocked. If a white pellet is visible, spin down the eluates and transfer the clear supernatant without a white pellet into a new reaction tube.
	<p><i>Buffer NTC</i></p> <ul style="list-style-type: none"> • Buffer NTI is used for double stranded DNA or long single stranded DNA above 500 bp in length. For shorter single stranded DNA or RNA, Buffer NTC is recommended. This buffer can be used for both, reaction clean up and gel extraction. • Buffer NTC can be ordered at MACHEREY-NAGEL (REF 740654.100)
Purification of RNA or single stranded DNA	

6.2 Ordering information

Product	REF	Pack of
NucleoSpin® Gel and PCR Clean-up	740609.10/.50/.250	10/50/250
Buffer NTI	740305.120	200 mL
Buffer NTB	740595.150	150 mL
Buffer NTC	740654.100	100 mL
Buffer NT3 (Concentrate) (for 125 mL Buffer NT3)	740598	25 mL
Collection Tubes (2 mL)	740600	1000
NucleoVac 24 Vacuum Manifold	740299	1
NucleoVac Mini Adapter	740297.100	100
NucleoVac Valves	740298.24	24
NucleoTrap®	740584.10/.50/.250	10/50/250
NucleoTraP®CR	740587.10/.50/.250	10/50/250

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

6.3 References

Vogelstein B., and D. Gillespie. 1979. Preparative and analytical purification of DNA from agarose. *Proc. Natl. Acad. Sci. USA* 76: 615–619.

6.4 Product use restriction / warranty

NucleoSpin® Gel and PCR Clean-up XS 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

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

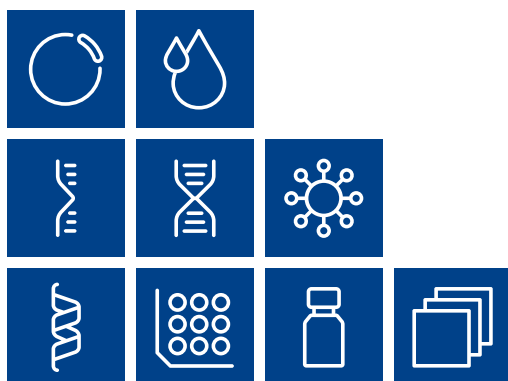
Trademarks:

NanoDrop® is a registered trademark of NanoDrop Technologies, Inc.

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

Roti® is a registered trademark of Carl Roth GmbH & Co. KG

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.



www.mn-net.com

MACHEREY-NAGEL



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

DE Tel.: +49 24 21 969-0
CH Tel.: +41 62 388 55 00
FR Tel.: +33 388 68 22 68
US Tel.: +1 484 821 0984

info@mn-net.com
sales-ch@mn-net.com
sales-fr@mn-net.com
sales-us@mn-net.com



A0xxxx/xxxx.xx