



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. ©2019

Genomic DNA from plant

User manual

NucleoMag[®] 384 Plant

June 2018 / Rev. 01

Contact MN

Germany and international

MACHEREY-NAGEL GmbH & Co. KG

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

Tel.: +49 24 21 969-0

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

Fax: +49 24 21 969-199

E-mail: info@mn-net.com

Technical Support Bioanalysis

Tel.: +49 24 21 969-270

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

USA

MACHEREY-NAGEL Inc.

2850 Emrick Blvd. · Bethlehem, PA 18020 · USA

Tel.: +1 484 821 0984

Toll-free: 888 321 6224 (MACH)

Fax: +1 484 821 1272

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

France

MACHEREY-NAGEL SARL à associé unique

1, rue Gutenberg · 67722 Hoerdt · France

Tel.: +33 388 68 22 68

Fax: +33 388 51 76 88

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

Switzerland

MACHEREY-NAGEL AG

Hirsackerstr. 7 · 4702 Oensingen · Switzerland

Tel.: +41 62 388 55 00

Fax: +41 62 388 55 05

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

www.mn-net.com

Table of contents

1 Components	4
1.1 Kit contents	4
1.2 Reagents, consumables, and equipment to be supplied by user	5
1.3 About this user manual	5
2 Product description	6
2.1 The basic principle	6
2.2 Kit specifications	6
2.3 Magnetic separation systems	6
2.4 Adjusting the shaker settings	7
2.5 Handling of beads	7
2.6 Storage and homogenization of samples	8
2.7 Elution procedures	9
3 Storage conditions and preparation of working solutions	9
4 Safety instructions	10
5 Protocol for the isolation of genomic DNA from plant samples in a 384-well format	12
5.1 Protocol at a glance	12
5.2 Detailed protocol	15
6 Appendix	18
6.1 Troubleshooting	18
6.2 Ordering information	19
6.3 Product use restriction / warranty	20

1 Components

1.1 Kit contents

NucleoMag[®] 384 Plant		
REF	1x 384 preps 744402.1	4 x 384 preps 744402.4
NucleoMag [®] V-Beads	2 x 1.5 mL	5 x 1.5 mL
Lysis Buffer MC1	2 x 60 mL	500 mL
Binding Buffer MC2	50 mL	80 mL
Wash Buffer MC3	75 mL	300 mL
Wash Buffer MC4	75 mL	300 mL
Wash Buffer MC5	60 mL	250mL
Elution Buffer MC6	30 mL	125 mL
RNase A (lyophilized)*	2 x 30 mg	7 x 30 mg
User manual	1	1

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

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

Reagents

80 % ethanol

Equipment / Consumables

Product	REF	Pack of
Lysis tubes for incubation of samples and lysis, e.g., Rack of Tubes Strips (1 set consists of 1 Rack, 12 Strips with 8 tubes (1.2 mL wells) each, and 12 Cap Strips)	740477 740477.24	4 sets 24 sets
Lysis plate for incubation of samples and lysis, e.g., Round well block (1 set consists of 1 Plate (96 x 1.2 mL round wells) each, and 12 Cap Strips)	740475 740475.24	4 sets 24 sets
Bead Tubes G for mechanical sample disruption, e.g., NucleoSpin® Bead Tubes G (50 tubes)	740817.50	50
Elution plate for collecting purified nucleic acids, e.g., Elution Plate U-bottom (96-well 0.3 mL microtiterplate with 300 µL u-bottom wells)	740486.24	24
e.g., Elution Plate Flat-bottom (96-well 0.3 mL microtiterplate with 300 µL flat-bottom wells)	740673	20

1.3 About this user manual

It is strongly recommended for first time users to read the detailed protocol sections of the **NucleoMag® 384 Plant** kit before using this product. Experienced users, however, may refer to the Protocol at a glance instead. The Protocol at a glance is designed to be used only as a supplemental tool for quick referencing while performing the purification procedure.

All technical literature is available online at www.mn-net.com.

2 Product description

2.1 The basic principle

The **NucleoMag[®] 384 Plant** procedure is based on reversible adsorption of nucleic acids to paramagnetic beads under appropriate buffer conditions. Plant tissue is extracted with CTAB-Lysis Buffer MC1. Adjusting the binding conditions of nucleic acid with Binding Buffer MC2 and addition of paramagnetic beads can be carried out simultaneously. After magnetic separation and removal of supernatant, the paramagnetic beads are washed with Wash Buffers MC3, MC4, and 80 % ethanol to remove contaminants and salt. There is no need for a drying step as ethanol from previous wash steps is removed by Wash Buffer MC5. Finally, highly purified DNA is eluted with low-salt Elution Buffer MC6 and can directly be used for downstream applications. The **NucleoMag[®] 384 Plant** is designed for automated high-throughput processing on standard liquid handling instruments in a 384-well format.

2.2 Kit specifications

NucleoMag[®] 384 Plant is designed for an automated preparation of DNA from plant samples in a 384-well format. The kit is designed for use with static magnetic separator plates. The purified DNA can be used directly as template for qPCR, next generation sequencing, or any kind of enzymatic reactions.

NucleoMag[®] 384 Plant allows easy automation on common liquid handling instruments. The actual processing time depends on the configuration of the instrument and the magnetic separation system used.

2.3 Magnetic separation systems

For the **NucleoMag[®] 384 Plant** kit, the use of a static magnetic separator plate in combination with a 96- or 384-multichannel device is recommended. Separation is carried out in a 384-well plate (not provided in the kit). The kit is also suitable for common 384-well plate magnetic separators

Magnetic separator	Separation plate or tube
Magnetic separator VP 771G-4A (V&P Scientific, Inc.)	384-well plates (deep well), (Greiner BioOne)
Alpaqua 384 Post Magnet Plate (Alpaqua Engineering, LLC)	

Static magnetic pins

Separators with static magnetic pins are recommended in combination with a suitable microplate shaker for optimal resuspension of the beads during the washing and elution steps. Alternatively, beads can be resuspended in the buffer by pipetting up and down several times. For fully-automated use on liquid handling workstations, a gripper tool is required, the plate is transferred to the magnetic separator for separation of the beads and transferred to the shaker module for resuspension of the beads.

2.4 Adjusting the shaker settings

When using a plate shaker for the washing and elution steps, the speed settings have to be adjusted carefully for each specific separation plate and shaker to prevent cross-contamination from well to well. Proceed as follows:

Adjusting shaker speed for binding and wash steps:

- Load 105 μL (for checking the binding step settings) or 80 μL (for checking the washing steps settings) dyed water to the wells of the separation plate. Place the plate on the shaker and start shaking with a moderate speed setting for 30 seconds. Turn off the shaker and check the plate surface for small droplets of dyed water.
- Increase speed setting, shake for an additional 30 seconds, and check the plate surface for droplets again.
- Continue increasing the speed setting until you observe droplets on top of the separation plate. Reduce speed setting, check again, and use this setting for the washing step.

Adjusting shaker speed for the elution step:

Load 100 μL dyed water to the wells of the collection plate and proceed as described above.

2.5 Handling of beads

Distribution of beads

A homogeneous distribution of the magnetic beads to the individual wells of the separation plate is essential for a high well-to-well consistency. Therefore, before distributing the beads, make sure that the beads are completely resuspended. Shake the storage bottle well or place it on a vortexer shortly. Premixing magnetic beads with the binding buffer allows easier homogenous distribution of the beads to the individual wells of the separation plate. During automation, a premix step before aspirating the beads/binding buffer mixture from the reservoir is recommended to keep the beads resuspended.

Magnetic separation time*

Attraction of the magnetic beads to the magnetic pins depends on the magnetic strength of the magnetic pins, the selected separation plate, distance of the separation plate from the magnetic pins, and the volume to be processed. The individual times for complete attraction of the beads to the magnetic pins should be checked and adjusted on each system. It is recommended using the separation plates or tubes specified by the supplier of the magnetic separator.

* Contact MN Technical Service for optimized program files and support protocols.

Washing the beads

Washing the beads can be achieved by shaking or mixing. In contrast to mixing by pipetting up and down, mixing by shaker or magnetic mixing allows simultaneous mixing of all samples. This reduces the time and number of tips needed for the preparation. Resuspension by pipetting up and down, however, is more efficient than mixing by a shaker or magnetic mix.

Method	Resuspension efficiency	Speed	Number of tips needed
Shaker	++	++	Low
Pipetting	+++	+*	High

+: acceptable, ++: good, +++: excellent

* 8-channel pipetting device

2.6 Storage and homogenization of samples

We recommend the use of young plant samples and to keep plants for about 12 h in the dark before collecting samples in order to reduce polysaccharide content.

Plant samples can be stored frozen, under ethanol or lyophilized. In many cases lyophilized, dried material can be easier processed and gives higher yield. If using dried samples, reduce the amount of starting material by the factor 5 (e.g., use 10 mg dried plant leaves instead of 50 mg fresh weight).

As plant tissue is very robust, the lysis procedure is most effective with well homogenized, powdered samples. Suitable methods include grinding with pestle and mortar in the presence of liquid nitrogen or using steel beads. We also recommend the use of other commercial homogenizers, bead mills, for example, Crush Express for 96-well homogenization (contact Saaten-Union Resistenzlabor GmbH, D-33818 Leopoldshohe), Tissue Striker (www.KisanBiotech.com) or Geno/Grinder 2000 (www.spexcsp.com or for Germany www.c3-analysentechnik.de).

Methods to homogenize samples

Homogenize samples in NucleoSpin® Bead Tubes G, Rack of Tube Strips or a Round well block with Cap Strips (see ordering Informations 6.2) filled with e.g. steel beads.

Alternatively homogenizing of larger samples can be performed by VA steel beads (diameter: 7 mm): Put 4–5 beads and plant material together into a 15 mL plastic tube (Falcon), chill the tube in liquid nitrogen and vortex for about 30 seconds (e.g., with a Multi Pulse Vortexer, contact Schütt Labortechnik GmbH, Postfach 3454, D-37024 Göttingen, Germany). Repeat this chilling and vortexing procedure until the entire plant material is ground to a powder. Chill the tube once more and remove the beads by rolling them out gently or with a magnet. Keep the material frozen throughout the whole homogenization procedure. Do not add nitrogen to the tube! This leads to sticking and loss of plant material attached to the beads.

2.7 Elution procedures

Purified DNA can be eluted directly with the supplied Elution Buffer MC6. Elution can be carried out in a volume of $\geq 50 \mu\text{L}$. It is essential to cover the NucleoMag[®] Beads completely with elution buffer during the elution step. The volume of dispensed elution buffer depends on the magnetic separation system (e.g., the position of the pellet inside the separation plate). For efficient elution, the magnetic bead pellet should be resuspended completely in the elution buffer. For some separators, higher elution volumes might be necessary to cover the whole pellet.

Elution is possible at room temperature. Yield can be increased by 15–20% if elution is performed at 55 °C.

3 Storage conditions and preparation of working solutions

Attention: Buffers MC3 and MC4 contain chaotropic salt! Buffer MC2 is highly flammable and irritant. Wear gloves and goggles!

Storage conditions:

- All components of the **NucleoMag[®] 384 Plant** kit should be stored at room temperature (18–25 °C) and are stable for up to one year.
- All buffers are delivered ready-to-use.

Before starting any **NucleoMag[®] 384 Plant** protocol, prepare the following:

- **RNase A:** Before first use, add the indicated volume of water to each vial of the lyophilized RNase A. **Store RNase A at 4 °C.**
- **80% ethanol:** Use molecular biology grade ethanol, dilute with appropriate water to 80%.

NucleoMag [®] 384 Plant		
REF	1 x 384 preps 744402.1	4 x 384 preps 744402.4
RNase A (lyophilized)	2 x 30 mg Add 2.5 mL water to each vial	7 x 30 mg Add 2.5 mL water to each vial





4 Safety instructions

The following components of the **NucleoMag® 384 Plant** kits contain hazardous contents.

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

GHS classification

Only harmful features do not need to be labeled with H and P phrases up to 125 mL or 125 g. *Mindergefährliche Eigenschaften müssen bis 125 mL oder 125 g nicht mit H- und P-Sätzen gekennzeichnet werden.*

Component	Hazard contents	GHS symbol	Hazard phrases	Precaution phrases
<i>Inhalt</i>	<i>Gefahrstoff</i>	<i>GHS-Symbol</i>	<i>H-Sätze</i>	<i>P-Sätze</i>
MC2	2-propanol 95–100 % <i>2-propanol 95–100 %</i> CAS 67-63-0	 DANGER GEFAHR	225, 319, 336	210, 233, 260D, 280sh
MC3	Ethanol 20–35 % <i>Ethanol 20–35 %</i> CAS 64-17-5	 WARNING ACHTUNG	226	210
MC4	Ethanol 20–35 % <i>Ethanol 20–35 %</i> CAS 64-17-5	 WARNING ACHTUNG	226	210
RNase A	RNase A, lyophilized 90–100 % <i>RNase, lyophilisiert 90–100 %</i> CAS 9001-99-4	 DANGER GEFAHR	317, 334	261sh, 342+311

The symbol shown on labels refers to further safety information in this section.

Das auf Etiketten dargestellte Symbol weist auf weitere Sicherheitsinformationen dieses Kapitels hin.

For further information please see Material Safety Data Sheets (www.mn-net.com).

Weiterführende Informationen finden Sie in den Sicherheitsdatenblättern (www.mn-net.com).

Hazard phrases

- H 225 Highly flammable liquid and vapour.
Flüssigkeit und Dampf leicht entzündbar.
- H 226 Flammable liquid and vapour.
Flüssigkeit und Dampf entzündbar.
- H 317 May cause an allergic skin reaction.
Kann allergische Hautreaktionen verursachen.
- H 319 Causes serious eye irritation.
Verursacht schwere Augenreizung.

- H 334 May cause allergy or asthma symptoms or breathing difficulties if inhaled.
Kann bei Einatmen Allergie, asthmaartige Symptome oder Atembeschwerden verursachen.
- H 336 May cause drowsiness or dizziness.
Kann Schläfrigkeit und Benommenheit verursachen.

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 260D Do not breathe vapors.
Dampf nicht einatmen.
- P 261sh Avoid breathing dust / vapors.
Einatmen von Staub / Dampf vermeiden.
- P 280sh Wear protective gloves / eye protection / face protection.
Schutzhandschuhe / Augenschutz tragen.
- P 342+311 If experiencing respiratory symptoms: Call a POISON CENTER / doctor / ...
Bei Symptomen der Atemwege: GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.

5 Protocol for the isolation of genomic DNA from plant samples in a 384-well format

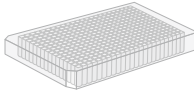
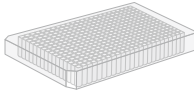
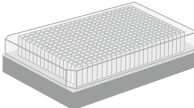
5.1 Protocol at a glance

For additional equipment and hardware requirements, refer to section 1.2 and 2.3 respectively.

For detailed information on each step, see page 15.

Before starting the preparation:

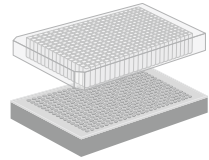
Check if RNase A was prepared according to section 3.

1 Homogenize and lyse plant sample material (~30 mg)	Add 200 μL MC1 and 10 μL RNase A Mix 56 °C, 60 min	
2 Clear lysates by centrifugation, transfer 50 μ L of cleared lysate for further processing	Centrifuge at 5,600 x g, 3 min Transfer 50 μL cleared lysate	
3 Bind DNA to NucleoMag[®] V-Beads	Add 4 μL NucleoMag[®] V-Beads and 50 μL MC2 <i>(Note: Resuspend the NucleoMag[®] V-Beads before removing them from the storage bottle)</i>	
	Mix by shaking for 5 min at RT <i>(Optional: Mix by pipetting up and down)</i>	
	Remove supernatant after 3 min separation	

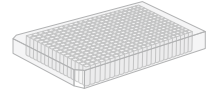
4 Wash with MC3

Remove separation plate from magnetic separator

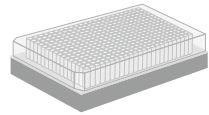
Add 80 µL MC3



Resuspend: Shake 5 min at RT
(Optional: Mix by pipetting up and down)



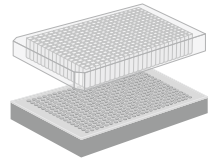
Remove supernatant after 2 min separation



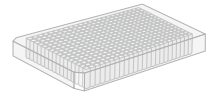
5 Wash with MC4

Remove separation plate from magnetic separator

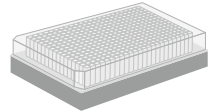
Add 80 µL MC4



Resuspend: Shake 5 min at RT
(Optional: Mix by pipetting up and down)



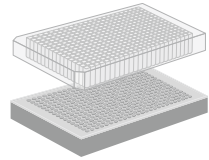
Remove supernatant after 2 min separation



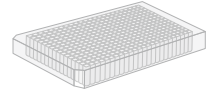
6 Wash with 80 % ethanol

Remove separation plate from magnetic separator

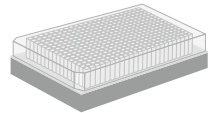
Add 80 μ L 80 % ethanol



Resuspend: Shake 5 min at RT
(Optional: Mix by pipetting up and down)



Remove supernatant after 2 min separation



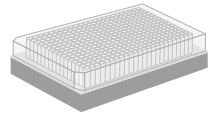
7 Wash with MC5

Leave Square-well Block on the magnetic separator!

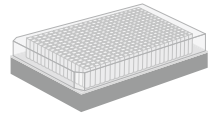
Add 80 μ L MC5

Incubate for 45–60 s

Note: Do not resuspend the beads in Buffer MC5!



Remove supernatant

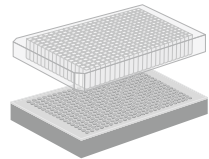


8 Elute DNA

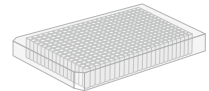
Remove separation plate from magnetic separator

Add 60 μ L MC6

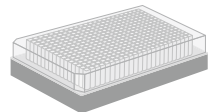
(Elution volume possible from 50 - 200 μ L; Optional: Elute at 55 °C)



Shake 10 min at RT
(Optional: Mix by pipetting up and down)



Separate 2 min and transfer DNA (supernatant) into elution plate/ tubes



5.2 Detailed protocol

This protocol is designed for magnetic separators with static pins and suitable plate shakers (see section 2.3). It is recommended to use a suitable 384-well separation plate (see section 2.3). Alternatively, isolation of DNA can be performed in other formats (see section 6.2). This protocol is for manual use and serves as a guideline for adapting the kit to robotic instruments.

Before starting the preparation:

Check if RNase A was prepared according to section 3.

1 Homogenize and lyse sample material

Homogenize about **30 mg (fresh)** or **< 10 mg (lyophilized) plant tissue**, for example, using microtube strips in a mixer mill, and add **200 µL Buffer MC1**. Do not moisten the rim. Close the individual wells with cap strips. Mix by vigorous shaking for 15–30 s. Spin briefly for 30 s at 1,500 x g to collect any sample from the cap strips. Incubate the closed strips at **56 °C for 60 min**.

Optional: If samples contain large amounts of RNA, we recommend the addition of 10 µL RNase A solution (stock solution 12 mg/mL) to the MC1 lysis mixture.

2 Clear lysates

Centrifuge the samples for **3 min** at a full speed (**5,600–6,000 x g**). Remove cap strips.

Transfer **50 µL of the cleared lysate** (equilibrated to room temperature) to a suitable separation plate. Do not moisten the rims of the well.

Note: See recommendations for suitable plates or tubes and compatible magnetic separators section 1.2.

3 Bind DNA to NucleoMag® V-Beads

Add **4 µL of NucleoMag® V-Beads** and **50 µL Buffer MC2** to each well of the separation plate. Mix by pipetting up and down 6 times and **shake for 5 min** at room temperature. Alternatively, when processing the kit without a shaker, pipette up and down 10 times and incubate for 5 min at room temperature.

Note: Be sure to resuspend the NucleoMag® V-Beads before removing them from the storage bottle. Vortex storage bottle briefly until a homogenous suspension has been formed.

Separate the magnetic beads against the side of the wells by placing the separation plate on a suitable magnetic separator. Wait at least **3 min** until all the beads have been attracted to the magnets. Remove and discard supernatant by pipetting.

Note: Do not disturb the attracted beads while aspirating the supernatant. The magnetic pellet is not visible in this step. Remove supernatant from the opposite side of the well.

4 Wash with MC3

Remove the separation plate from the magnetic separator.

Add **80 µL Buffer MC3** to each well and resuspend the beads by shaking until the beads are resuspended completely (**5 min**). Alternatively, resuspend beads completely by repeated pipetting up and down (15 times).

Separate the magnetic beads by placing the separation plate on the magnetic separator. Wait at least **2 min** until all the beads have been attracted to the magnet. Remove and discard supernatant by pipetting.

5 Wash with MC4

Remove the separation plate from the magnetic separator.

Add **80 µL Buffer MC4** to each well and resuspend the beads by shaking until the beads are resuspended completely (**5 min**). Alternatively, resuspend beads completely by repeated pipetting up and down (15 times).

Separate the magnetic beads by placing the separation plate on the magnetic separator. Wait at least **2 min** until all the beads have been attracted to the magnet. Remove and discard supernatant by pipetting.

6 Wash with 80 % ethanol

Remove the separation plate from the magnetic separator.

Add **80 µL 80 % Ethanol** to each well and resuspend the beads by shaking until the beads are resuspended completely (**5 min**). Alternatively, resuspend beads completely by repeated pipetting up and down (15 times).

Separate the magnetic beads by placing the separation plate on the magnetic separator. Wait at least **2 min** until all the beads have been attracted to the magnet. Remove and discard supernatant by pipetting.

7 Wash with MC5

Leave the separation plate on the magnetic separator.

Note: Supernatant is colorless, magnetic bead pellet is clearly visible.

Gently add **80 µL Buffer MC5** to each well and incubate for **45–60 s** while the beads are still attracted to magnets. Then aspirate and discard the supernatant.

Note: Do not resuspend the beads in Wash Buffer MC5. This step is to remove traces of ethanol and eliminates a drying step!

8 Elution

Remove the separation plate from the magnetic separator.

Add desired volume of **Buffer MC6 (50–200 µL)** to each well of the separation plate and resuspend the beads by shaking **5–10 min** at room temperature. Alternatively, resuspend beads completely by repeated pipetting up and down and incubate for **5–10 min** at room temperature.

Separate the magnetic beads by placing the separation plate on the magnetic separator. Wait at least **2 min** until all the beads have been attracted to the magnets. Transfer the supernatant containing the purified genomic DNA to the Elution Plate.

Note: Yield can be increased by 15–20 % by using pre-warmed elution buffer (55 °C) or by incubating the bead / elution buffer suspension at 55 °C for 10 min.

6 Appendix

6.1 Troubleshooting

Problem	Possible cause and suggestions
Poor DNA yield	<i>Elution buffer volume insufficient</i>
	Beads pellet must be covered completely with elution buffer
	<i>Insufficient performance of elution buffer during elution step</i>
	Remove residual buffers during the separation steps completely. Remaining buffers decrease efficiency of following wash steps and elution step.
	<i>Beads dried out</i>
Do not let the beads dry as this might result in lower elution efficiencies.	
<i>Partial elution in Wash Buffer MC5 already</i>	
Keep the beads on the magnet while dispensing Wash Buffer MC5. Do not resuspend beads in this buffer, and do not incubate beads in this buffer for more than 2 min, as this buffer is water-based and might elute the DNA already.	
<i>Aspiration of attracted bead pellet</i>	
Do not disturb the attracted beads while aspirating the supernatant, especially when the magnetic pellet is not visible in the lysate.	
<i>Incubation after dispensing beads to lysate</i>	
Mix immediately after dispensing NucleoMag® V-Beads/Buffer MC2 to the lysate.	
Low purity	<i>Insufficient washing procedure</i>
	Use only the appropriate combinations of separator and plate, for example.
Suboptimal performance of DNA in downstream applications	<i>Carry-over of ethanol from 80 % ethanol wash solution</i>
	Be sure to remove all of the ethanolic wash solution, as residual ethanol interferes with downstream applications.
	<i>Low purity</i>
See above	

Problem	Possible cause and suggestions
Carry-over of beads	<i>Time for magnetic separation too short</i>
	Increase separation time to allow the beads to be completely attracted to the magnetic pins before aspirating any liquid from the well.
	<i>Aspiration speed too high (elution step)</i>
	High aspiration speed during the elution step may cause bead carry-over. Reduce aspiration speed for elution step.
Cross contamination	<i>Contamination of the rims</i>
	Do not moisten the rims of the processing plate when transferring the plant lysate. If the rim of the wells is contaminated, seal the processing plate with Self-adhering PE Foil (see ordering information) before starting the shaker.

6.2 Ordering information

Product	REF	Pack of
NucleoMag [®] 384 Plant	744402.1	1 x 384 preps
	744402.4	4 x 384 preps
NucleoMag [®] SEP Just for use with 96-well plates, such as 740481 (Deep-well plate). Not recommended for the use of 384-well blocks.	744900	1
Square-well Blocks (96-well blocks with square wells for use with NucleoMag SEP 744900)	740481	4
	740481.24	24
Self-adhering PE Foil	740676	50 sheets
Rack of Tube Strips (set consists of 1 Rack, 12 Tube Strips with 8 tubes each, and 12 Cap Strips)	740477	4 sets
	740477.24	24 sets
Round well block with Cap Strips (set consists of 1 Plate (96 x 1.2 mL round wells) each, and 12 Cap Strips)	740475	4 sets
	740475.24	24 sets
NucleoSpin [®] Bead Tubes G (2 mL tubes with steel beads for sample homogenization)	740817.50	50
Elution Plate U-bottom	740486.24	24
Elution Plate Flat-bottom	740673	20
Cap Strips	740638	30 strips

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

6.3 Product use restriction/warranty

NucleoMag® 384 Plant 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

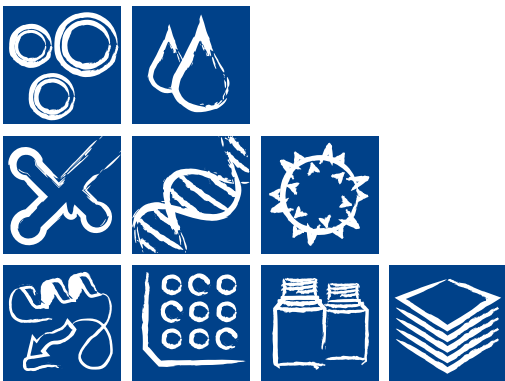
Trademarks:

NucleoMag is a registered trademark of MACHEREY-NAGEL 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.



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



www.mn-net.com

MACHEREY-NAGEL



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

DE / International:

Tel.: +49 24 21 969-0
 Fax: +49 24 21 969-199
 E-mail: info@mn-net.com

CH:

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

FR:

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

US:

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

