

NucleoMag® RNA/ DNA Pro – MagnetaPure 32+

Protocol details

Application	Isolation of RNA and DNA from tissue and plant leaf material
Kit	NucleoMag® RNA Pro
REF	744370.1
Protocol name	NMRNADNAPROaR1, NMRNADNAPRObR1



Eight easy steps

Procedure	
1	Perform lysis according to the user manual NucleoMag® RNA Pro.
2	Fill the 96-well Deep-well plates 1 & 2 according to the tables for DNA and RNA sections below.
3	Load the plates on the MagnetaPure 32+.
4	Insert tip combs on the mounting grooves. *
5	Select the protocol from the instrument menu and start protocol NMRNADNAPROaR1.
6	Remove the plates after the DNA extraction is finished.
7	For rebinding transfer 300 µL of the rDNase reaction mixture containing the RNA from plate 1 column 6 + 12 to the second plate (column 1 + 7), Please refer to the loading table and schema Plate 2 (RNA section)**
8	Place Plates 2 into the MagnetaPure 32+ and proceed immediately with the RNA extraction and proceed with protocol NMRNADNAPRObR1.

Note: * Please equip all tip combs in order to cover the magnetic rods in used and unused wells.

** To remove the plate the tip combs have to be temporarily removed, they have to be reused in their respective positions, please ensure that the attached beadpellet is not disturbed during this..

Loading table for plate (DNA Section)

Position	Reagents	Samples per plate
Column 1 + 7	Cleared Lysate (350 µL), Binding Reagent (250 µL)*, NucleoMag® B-Beads (20 µL)	Sample 1-8 Sample 9-16
Column 2 + 8	Wash Buffer MRW (900 µL)	Sample 1-8 Sample 9-16
Column 3 + 9	Wash Buffer DNA Wash (900 µL)	Sample 1-8 Sample 9-16
Column 4 + 10	Elution Buffer DNA Elute (100 µl)	Sample 1-8 Sample 9-16
Column 5 + 11	Empty	Sample 1-8 Sample 9-16
Column 6 + 12	rDNase reaction mixture (300 µl)	Sample 1-8 Sample 9-16

Note: Please refer to the image below for a visual representation of the loading scheme

* Use binding reagent for different sample types according to the user manual NucleoMag® RNA Pro.



Loading scheme (DNA section)

	1	2	3	4	5	6	7	8	9	10	11	12	
A													
B													
C													
D													
E													
F													
G													
H													
	Binding (620 µL)	1 st Wash (900 µL)	2 nd Wash (900 µL)	DNA Elution (100 µL)		DNA digest (300 µL)	Binding (620 µL)	1 st Wash (900 µL)	2 nd Wash (900 µL)	DNA Elution (100 µL)		DNA digest (300 µL)	

Loading table for plate (RNA Section)

Position	Reagents	Samples per plate
Column 1 + 7	Rebinding Buffer MRB (350 µl) + rDNase reaction mixture containing RNA (300 µl)*	Sample 1-8 Sample 9-16
Column 2 + 8	Ethanol 70% (900 µL)	Sample 1-8 Sample 9-16
Column 3 + 9	Empty	Sample 1-8 Sample 9-16
Column 4 + 10	Empty	Sample 1-8 Sample 9-16
Column 5 + 11	Empty	Sample 1-8 Sample 9-16
Column 6 + 12	Elution Buffer MRE (100 µL)	Sample 1-8 Sample 9-16

Note: Please refer to the image below for a visual representation of the loading scheme

* The rDNase reaction mixture is transferred from column 6 & 12 of plate 1 after running the DNA section of the protocol. Plate 1 can be removed when the Instrument finished part a. Perform the transfer, place plate 2 into the instrument and continue with the protocol part b.

Loading scheme (RNA section)

	1	2	3	4	5	6	7	8	9	10	11	12	
A	Rebinding (350 µL + 300 µL)	3 rd Wash (900 µL)				RNA Elution (100 µL)		Rebinding (350 µL + 300 µL)				RNA Elution (100 µL)	
B													
C													
D													
E													
F													
G													
H													

Additional consumables and instrumentation

Product	Specification	REF
MagnetaPure 32+	Automated nucleic extraction system for MACHEREY-NAGEL's NucleoMag® kits enabling parallel processing of up to 32 samples	747010
96 Deep-well Plate	96 deep-well plates for MagnetaPure 32+ (25 pieces)	744955
Tip Combs	8-well tip combs for MagnetaPure 32+ (50 pieces)	744960

Disclaimer

Information

MACHEREY-NAGEL GmbH & Co. KG makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. MACHEREY-NAGEL GmbH & Co. KG cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your MACHEREY-NAGEL representative. This publication may contain reference to applications and products which are not available in all markets. Please check with your local sales representative. All mentioned trademarks are protected by law. 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). Reference herein to any specific commercial product, process, or service by trade name, trademark, manufacturer, or otherwise, does not necessarily constitute or imply its endorsement, recommendation, or support by MACHEREY-NAGEL GmbH & Co. KG. Any views or opinions expressed herein by the authors' do not necessarily state or reflect those of MACHEREY-NAGEL. NucleoMag® is a registered trademark of MACHEREY-NAGEL GmbH & Co. KG, Düren, Germany.