

NucleoMag[®] RNA Pro – MagnetaPure 32+

Protocol details

| Application | Highly pure RNA from tissue and plant leaf material |
|---------------|---|
| Kit | NucleoMag [®] RNA Pro |
| REF | 744360 (.1 /.4) |
| Protocol name | NMRNAPROaRev01, NMRNAPRObRev01 |



Seven easy steps

| Procedure | |
|-----------|---|
| 1 | Perform lysis according to the user manual NucleoMag [®] RNA Pro. |
| 2 | Fill the 96-well Deep-well plates according to the table 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 NMRNAPROaRev01. |
| 6 | After the run (approx. 44 min), dispense 350 µL of Binding Buffer MRB to the DNA digestion (column 4 + 10). ** |
| 7 | Place the plate back into the MagnetaPure 32+ slide in the tip combs and proceed with protocol NMRNAPRObRev01 (approx. 34 min). |

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 can be reused in their respective positions, but we recommend using a new set of tip combs after placing the plate back into the instrument.

Loading table

| 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 | Ethanol 70% (900 µL) | Sample 1-8 Sample 9-16 |
| Column 4 + 10 | rDNase reaction mixture (300 µL) | Sample 1-8 Sample 9-16 |
| Column 5 + 11 | Ethanol 70% (900 µL) | 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

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

Loading scheme

| | 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 digest (300 µL + 350 µL) | 3 rd Wash (900 µL) | Elution (100 µL) | Binding (620 µL) | 1 st Wash (900 µL) | 2 nd Wash (900 µL) | DNA digest (300 µL + 350 µL) | 3 rd Wash (900 µL) | Elution (100 µL) | |

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.