



NucleoSpin® Gel and PCR Clean-up

Kit contents

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. Clontech and the MN logo are trademarks and the property of Clontech Laboratories, Inc. ©2015

NucleoSpin® Gel and PCR Clean-up

| REF | 12 preps 740609.12C | 240 preps 740609.240C |
|---|---|---|
| Binding Buffer NT1 | 40 mL | 200 mL |
| Wash Buffer NT3 (Concentrate) | 6 mL (add 24 mL ethanol before first use) | 60 mL (add 240 mL ethanol before first use) |
| Elution Buffer NE | 30 mL | 125 mL |
| NucleoSpin® Gel and PCR Clean-up Columns (yellow rings) | 12 | 4 x 60 |
| User manual | 1 | 1 |

How to use the kit

Please see the protocol information how to use the kit (see next pages). For further questions and more detailed information, please contact MACHEREY-NAGEL at tech-bio@mn-net.com for protocol information how to use the kit on specific robotic instruments.

For storage conditions, product use restrictions, and safety information, please see the general NucleoSpin® Gel and PCR Clean-up user manual.

General information

| | |
|----------------------|--|
| Application: | Clean-up |
| Kit: | NucleoSpin® Gel and PCR Clean-up (REF 740609.240C) instead of: QIAquick® PCR Purification Kit |
| Sample material: | Amplification reaction |
| Protocol name: | QIAquick® PCR Purification protocol (Check on QIAcube® web portal for most recent protocol versions.) |
| Editable parameters: | Elution volume: 30–100 µL; default 50 µL Fill-up volume: 0–96 µL; default 50 µL |

Using the kit

1. Fill the designated buffer bottles with the buffers according to the buffer table on page 3.
2. Samples: 5–100 µL PCR reaction
For PCR sample volume < 100 µL, the QIAcube® has to fill up the sample volume to 100 µL. Make sure to enter the volume of Elution Buffer NE that the QIAcube® should use to adjust the sample volume to 100 µL before starting the purification procedure.
3. Place Sample Tubes (2 mL Safe-Lock microcentrifuge tubes) into sample rack (shaker).
4. Insert disposable Filter Tips 1000 µL and 200 µL.
5. General equipment setup is shown below.

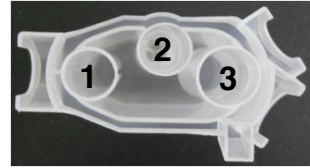
Additional materials

Refer to the QIAcube® protocol sheet for required consumables (e.g., sample tubes, collection tubes, instrument accessories, disposable tips, etc.) and software requirements.

Rotor adapter

| Position | Labware | Lid position |
|----------|---|--------------|
| 1 | NucleoSpin® Gel and PCR Clean-up Column | L1 |
| 2 | – | – |
| 3 | 1.5 mL collection tube* | L3 |

L1



L2

L3

* Sarstedt, Micro tube 1.5 mL Safety Cap

Buffers (Reagent Bottle Rack)

| Position | MN Reagent | Replaced QIAGEN® Reagent |
|----------|------------|--------------------------|
| 1 | Buffer NT1 | Buffer PB or Buffer PBI |
| 2 | – | – |
| 3 | – | – |
| 4 | – | – |
| 5 | Buffer NT3 | Buffer PE |
| 6 | Buffer NE | Buffer EB |

Note: MN is not addressing to use this kit on specific robots, it is customer's choice and interest how to use the kit for nucleic acid isolation.

NucleoSpin® kits on QIAcube®

The use of NucleoSpin® kits on the QIAcube® is at your own liability. MACHEREY-NAGEL is not responsible for loss of warranty claims or other consequences.

Trademarks: QIAGEN, QIAcube, and QIAamp are registered trademarks of the QIAGEN Group; NucleoSpin is a registered trademark of MACHEREY-NAGEL GmbH & Co. KG