



NucleoSpin® RNA

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. Unless otherwise indicated, all trademarks are the property of Clontech Laboratories, Inc. 80015

NucleoSpin® RNA		
REF	12 preps 740955.12c	240 preps 740955.240c
Lysis Buffer RA1	10 mL	250 mL
Buffer RA3 (Concentrate)	6 mL (add 24 mL ethanol before first use)	100 mL (add 400 mL ethanol before first use)
Membrane Desalting Buffer MDB	25 mL	250 mL
Reaction Buffer for rDNase	7 mL	30 mL
rDNase, RNase-free (lyophilized)	1 vial (size D) (add 130 µL RNase-free H ₂ O for reconstitution)	5 vials (size F) (add 550 µL RNase-free H ₂ O for reconstitution to each vial)
RNase-free H ₂ O	13 mL	125 mL
NucleoSpin® Filters (violet rings)	12	240
NucleoSpin® RNA Columns (light blue rings – plus Collection Tubes)	12	240
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 storage conditions, product use restrictions, and safety information, please see the general NucleoSpin® RNA user manual.

General information

Application:	RNA
Kit:	NucleoSpin® RNA (REF 740955.240c) instead of: RNeasy® Mini Kit
Sample material:	Animal cells cultured cells
Protocol name:	Purification of total RNA from animal cells (QIAshredder homogenization and on-column DNase digest) (Check on QIAcube® web portal for most recent protocol versions.)
Short protocol name:	QIAshredder DNase digest

Using the kit

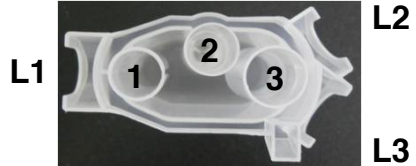
1. Fill the designated buffer bottles with the buffers according to buffer table on page 3.
2. Sample homogenization in a 2 mL Sample Tube.
Disrupt 5×10^6 – 1×10^7 cells in 600 μ L Buffer RA1 containing 6 μ L β -mercaptoethanol or 12 μ L of a 1 M TCEP stock solution (thaw cells before starting).
3. Place sample tubes into sample rack.
4. Fill required volume of rDNase reaction mix into a 2 mL Safe-Lock microcentrifuge tube.
5. Place rDNase reaction mix in position A of the Microcentrifuge Tube Slots.
6. Insert disposable Filter Tips 1000 μ L and 1000 μ L wide-bore.
7. General equipment setup is shown on page 3.

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® RNA Column	L1
2	NucleoSpin® Filters (no lid, violet ring)	–
3	1.5 mL collection tube*	L3



* Sarstedt, Micro tube 1.5 mL Safety Cap

Buffers (Reagent Bottle Rack)

Position	MN Reagent	Replaced QIAGEN® Reagent
1	–	–
2	70 % ethanol	70 % ethanol
3	–	–
4	Buffer MDB	Buffer RW1
5	Buffer RA3	Buffer RPE
6	RNase-free H ₂ O	RNase-free water

Microcentrifuge Tube Slots

	Position A	Position B	Position C
Content:	rDNase reaction mix	–	–
Tube:	2 mL Safe-Lock microcentrifuge tube	–	–

Required volume of rDNase reaction mix in Microcentrifuge Tube Slots

No. of samples	Diluted Carrier RNA (Microcentrifuge Tube Slot C)
2	213 µL (21 µL DNase + 192 µL DNase Reaction Buffer)
3	300 µL (32 µL DNase + 268 µL DNase Reaction Buffer)
4	386 µL (43 µL DNase + 343 µL DNase Reaction Buffer)
5	472 µL (54 µL DNase + 418 µL DNase Reaction Buffer)
6	559 µL (65 µL DNase + 494 µL DNase Reaction Buffer)
7	645 µL (76 µL DNase + 569 µL DNase Reaction Buffer)
8	731 µL (86 µL DNase + 645 µL DNase Reaction Buffer)
9	818 µL (97 µL DNase + 721 µL DNase Reaction Buffer)
10	904 µL (108 µL DNase + 796 µL DNase Reaction Buffer)
12	1077 µL (130 µL DNase + 947 µL DNase Reaction Buffer)

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