

# NucleoMag<sup>®</sup> SEP Magnetic Separator

## User manual

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](http://takarabio.com)

For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Takara Bio Inc. ©2018

April 2017 / Rev. 03

## Table of contents

1	Contents	4
2	Product description	4
3	Use of NucleoMag® SEP	4
3.1	Recommended separation plates	4
3.2	Separation of beads	5
3.3	NucleoMag® SEP care	5
4	Appendix	6
4.1	Ordering information	6
4.2	Product use restriction / warranty	6

# 1 Contents

NucleoMag® SEP	
REF	744900
NucleoMag® SEP Magnetic Separator	1
User manual	1

## 2 Product description

The NucleoMag® SEP Magnetic Separator is designed for the rapid manual purification of nucleic acids using the NucleoMag® 96 kits. The NucleoMag® kits are based on MACHEREY-NAGEL's magnetic-bead technology and allow rapid simultaneous purification of high-quality nucleic acids.

The NucleoMag® SEP can also be used on common liquid handling instruments, allowing a complete automation of the NucleoMag® kits in combination with a gripper tool.

## 3 Use of NucleoMag® SEP

### 3.1 Recommended separation plates

For optimal performance of the NucleoMag® SEP together with the NucleoMag® kits the following separation plates are recommended:


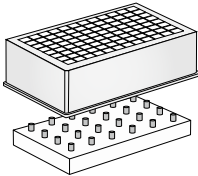
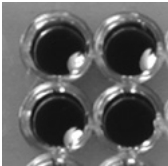
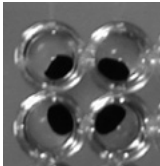
Elution Plate U-bottom: REF 740486.24

Square-well Block: REF 740481/.24

For more information refer to the appropriate kit protocol.

### 3.2 Separation of beads

During the washing and elution steps of the NucleoMag® procedure, the NucleoMag® Beads are freely dispersed in solution. Mixing of beads and buffer occurs by pipetting up and down or by shaking. To separate the beads from the solution, the separation plate is transferred onto the NucleoMag® SEP. The NucleoMag® Beads are attracted to the magnet, while the wash buffer can be removed from the well or the elution buffer containing the purified DNA can be recovered, respectively.

	During washing/elution step	At the end of washing/elution step
	Beads are mixed with buffer by pipetting up and down or by shaking.	Beads are attracted to the magnet so the wash buffer can be removed or the elution buffer with DNA can be recovered. Beads are held back in the well.
Position of separation plate	On work bench/robot deck or placed on shaker 	Transferred to NucleoMag® SEP 
NucleoMag® Beads	Freely dispersed in solution 	Attracted to magnet 

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 on each system. It is recommended to use the separation plates or tubes specified in the appropriate kit protocol.

### 3.3 NucleoMag® SEP care

The NucleoMag® SEP consists of a PTFE base and magnetic pins. Do not autoclave. If necessary, clean the NucleoMag® SEP by rinsing it with water or by wiping off dirt and liquids with a towel. NucleoMag® SEP is resistant to diluted acetic acid, ethanol, guanidine-HCl, NaCl, NaOH, SDS. However, prolonged exposure of the separator to these chemicals should be avoided.

## 4 Appendix

### 4.1 Ordering information

Product	REF	Pack of
NucleoMag® SEP Magnetic Separator (for use with 96-well plates)	744900	1
Square-well Block	740481	4
	740481.24	24
Elution Plate U-bottom	740486.24	24

Related products	REF	Pack of
NucleoMag® SEP Mini magnetic separator (for use with up to 12 x 1.5 mL or 2 mL reaction tubes)	744901	1
NucleoMag® SEP Maxi magnetic separator (for use with up to 4 x 50 mL tubes)	744902	1
NucleoMag® SEP 24 magnetic separator (for use with 24-well plates)	744903	1

### 4.2 Product use restriction / warranty

**NucleoMag® SEP** is 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

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



[www.mn-net.com](http://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](mailto:info@mn-net.com)

CH:

Tel.: +41 62 388 55 00  
 Fax: +41 62 388 55 05  
 E-mail: [sales-ch@mn-net.com](mailto: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](mailto:sales-fr@mn-net.com)

US:

Tel.: +1 484 821 0984  
 Fax: +1 484 821 1272  
 E-mail: [sales-us@mn-net.com](mailto:sales-us@mn-net.com)