



NucleoMag SEP Magnetic Separator

User manual

*Visit our website
for more details!
click here...*

This product is 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	PT5010-1 PR943198
For technical support 1.800.662.2566 tech@clontech.com	

February 2003/Rev. 01

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	4
3.3	NucleoMag SEP care	5
4	Ordering information	6
5	Product Use Restriction / Warranty	6

1 Contents

	NucleoMag SEP
Cat. No.	744 900
NucleoMag SEP Magnetic Separator	1
Protocol	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


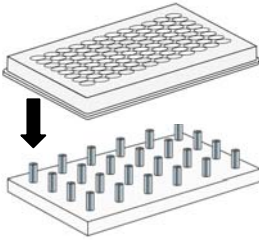
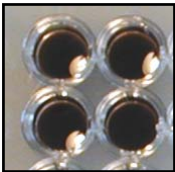

For optimal performance of the NucleoMag SEP together with the NucleoMag kits the following separation plates are recommended:

MN Elution Plate, U-Bottom	Cat. No. 744 672
MN Square-well Block	Cat. No. 744 670

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 teflon 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 Ordering information

Product	Cat. No.	Pack of
NucleoMag SEP Magnetic Separator	744 900	1
Square-well blocks	740 670	20
Elution Plate, U-bottom (to be used as separation plate)	740 672	20

5 Product Use Restriction / Warranty

NucleoMag SEP components were developed, designed and sold **for research purposes only**.

This MACHEREY-NAGEL product is shipped with documentation stating specifications and other technical information. MACHEREY-NAGEL guarantees 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 an extra copy.

MACHEREY-NAGEL does not warrant against 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; against defects in products or components not manufactured by MACHEREY-NAGEL, or against 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, PRODUCTIVITY, 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, incidental, 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.

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.

Please contact:

MACHEREY-NAGEL Germany
Tel.: +49-2421/969 270 and 275
e-mail: TECH-BIO@mn-net.com