

## NucleoVac 24 Vacuum Manifold

### 1 Contents

NucleoVac 24 Vacuum Manifold	
REF	740299
Vacuum manifold	1
NucleoVac Mini Adapter	24
Luer plugs	24
Tubing connection	2
Closing plug	1
Leaflet	1

### 2 Product description

The NucleoVac 24 Vacuum Manifold allows the parallel processing of silica columns for nucleic acid purification by vacuum. The silica columns (e.g., NucleoSpin® columns, NucleoSnap columns) can be attached to the upper luer slots of the vacuum manifold. Binding and washing steps of the NucleoSpin® or NucleoSnap procedures are carried out by vacuum instead of centrifugation, enabling time-saving procedures (NucleoSpin® products) and processing of large sample volumes (NucleoSnap products). Please see section 5 for an overview of NucleoSpin® kits suitable for vacuum processing. Support protocols are available online at [www.mn-net.com](http://www.mn-net.com). For processing NucleoSnap kits, please refer to the kit's user manual.

### 3 Storage conditions

All components should be stored at room temperature.

### 4 Safety instructions

When working with vacuum, it is recommended to wear protective goggles. For additional safety information on kit components, please refer to the kit's user manual.

### 5 Setup of the NucleoVac 24 Vacuum Manifold

See figures on reverse page

Close the lower threaded opening of the NucleoVac 24 Vacuum Manifold with the rubber closing plug ①. Screw in the tubing connection ② in the upper opening ③ manually. Do not use a wrench as the conical screw thread might crack if the connection is over-tightened. Attach tube (not included in the product).

*Note:* The product contains two different tubing connections for different tube diameters. Make sure to choose the tubing connection that enables complete tightness.

Connect NucleoVac 24 Vacuum Manifold to a vacuum source by attaching the tube to the vacuum source inlet.

*Note:* It is strongly recommended to use an overflow bottle between the NucleoVac 24 Vacuum Manifold and the vacuum source to avoid any contamination of the vacuum source, e.g., pump.

For processing NucleoSpin® ④ or NucleoSnap columns with the NucleoVac 24 Vacuum Manifold, insert NucleoSpin® or NucleoSnap columns into the Luer openings ⑤, a.

To avoid cross-contamination, it is recommended to use disposable NucleoVac Mini Adapters ⑥, see ordering information ⑥. For processing samples with different flow rates (e.g., large volume samples), it is recommended to use NucleoVac Stop-cocks ⑦, see ordering information ⑦) for maintaining constant vacuum pressure throughout the whole procedure ⑧. Close unused positions with Luer plugs ⑧, d.

### 6 Cleaning of the NucleoVac 24 Vacuum Manifold

Disconnect the NucleoVac 24 Vacuum Manifold and remove NucleoVac Mini Adapters and NucleoVac Stop-cocks. Discard NucleoVac Mini Adapters.

Remove the tubing connection and closing plug. Pour out liquid waste from the NucleoVac 24 Vacuum Manifold.

Rinse the NucleoVac 24 Vacuum Manifold with standard laboratory cleaning solutions and finally with water. If necessary, use standard disinfectant.

*Note:* The waste capacity of the manifold is 900 mL for non-foaming liquids and is reduced by foaming liquids. Empty the NucleoVac 24 manifold regularly!

Allow the NucleoVac 24 Vacuum Manifold and accessories to air-dry.

### 7 Ordering information

Product	REF	Pack of
NucleoVac 24 Vacuum Manifold	740299	1
NucleoVac Mini Adapter	740297.100	100
NucleoVac Stop-cock	740298.24	24

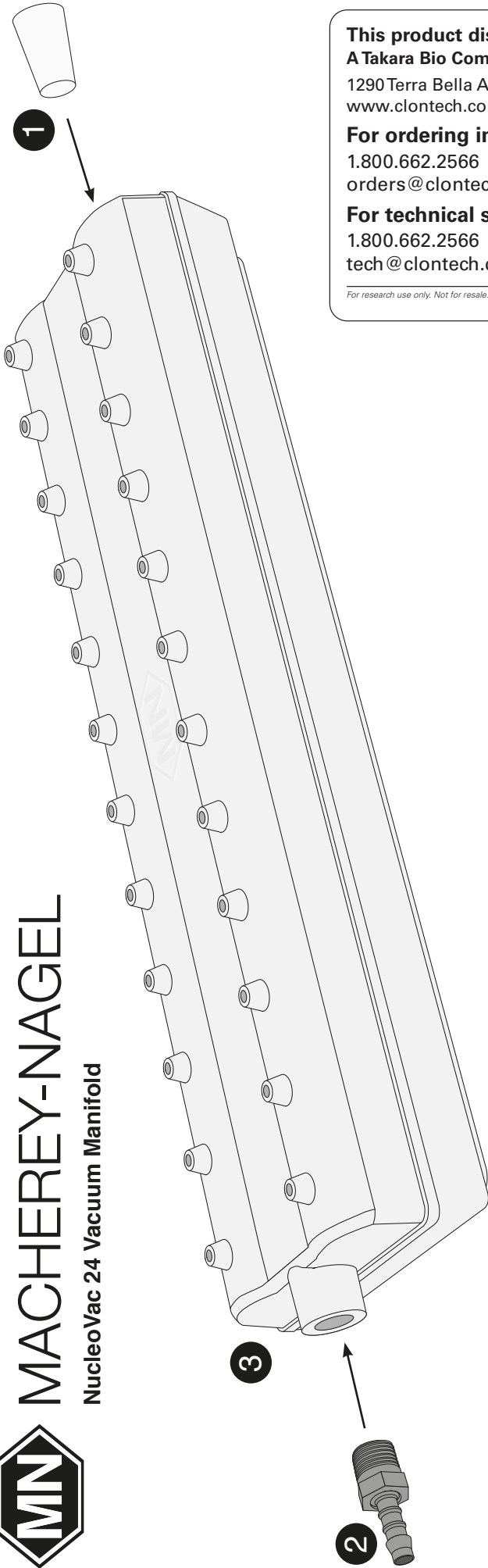
### 8 Product use restrictions / warranty

NucleoVac 24 Vacuum Manifold is developed, designed, and sold for research purposes only. We offer a one year warranty from date of delivery that our products will conform to applicable specifications set forth in the product specifications if not sold to persons set forth in § 13 of German Civil Code. In such a case, the provisions of the German Civil Code shall be valid. This warranty does not cover defects in or damage to the products which are due to improper installation or maintenance, misuse, neglect or any use other than ordinary commercial application. The seller shall not be liable for the goods being fit for a particular purpose unless otherwise agreed upon, to which the buyer intends to put them. This warranty is strictly exclusive. 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; REPRODUCIBILITY; DURABILITY; FITNESS FOR A PARTICULAR PURPOSE OR USE; MERCHANTABILITY; CONDITION; OR ANY OTHER MATTER WITH RESPECT TO MACHEREY-NAGEL PRODUCTS OR THE SALE OF MACHEREY-NAGEL PRODUCTS. If not superseded by these terms the current terms and conditions for the sale of goods (export version) are additionally valid.



# MACHEREY-NAGEL

NucleoVac 24 Vacuum Manifold



**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. ©2016

