

## Automated nucleic acid purification using MACHERY-NAGEL NucleoMag<sup>®</sup> kits

Freedom EVO<sup>®</sup> 75 optimized configuration

### Automation doesn't get easier than this!

If you have limited lab space and wish hand-free, fully automated magnetic-bead based purification of highly pure nucleic acid this affordable bench top workstation is the ideal solution.

### Ready-to-run optimized protocols

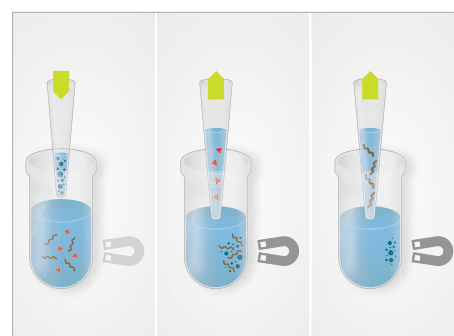
Optimized configurations and protocols have been developed in collaboration with MACHERY-NAGEL for a variety of sample types, including blood, tissue, bacteria and cells. The solution offers flexible processing of batches from 1 to 96 samples.

**Easy to use:** Nucleic Acid Purification dedicated TouchTools<sup>™</sup> interface

**Start instantly:** ready-to-run protocols included

**Superior quality:** proven yields and purity with MACHERY-NAGEL kits

**Convenient:** pre-configured system optimized to meet your needs



NucleoMag principle



Nucleic acid purification workstation

## Fast and efficient purification

Simplify your daily routine!

### What can I use it for?

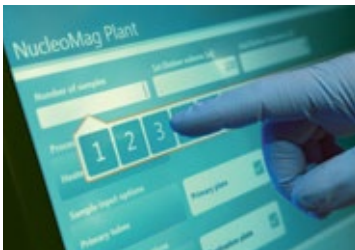
- Isolation of **genomic DNA from whole blood** with the NucleoMag Blood 200 µL kit
- Isolation of **genomic DNA from human or animal tissue, cultured cells, or bacteria** with the NucleoMag® 96 Tissue kit

### What is included?

- Two liquid handling channels
- Shaking and incubation module
- Support for plates and disposable tips
- Robotic manipulator for automatic sample transfer on and off the NucleoMag® SEP magnetic separator
- All necessary software and TouchTools interface

### Simple and intuitive

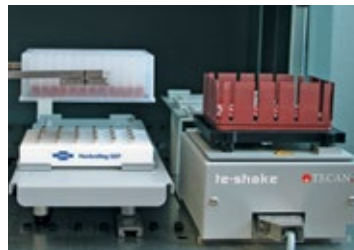
The TouchTools displays exactly what the operator needs for a run. User prompts with graphics provide step-by-step setup instructions for worktable set up, clearly and concisely.



Nucleid acid purification wizard

### Reliable optimized method

- In-well contained procedure minimizes risk of cross-contamination
- Small elution volumes ≥ 50 µL possible
- Yield does not depend on elution volume



Transport from magnet to shaker

### Sample traceability

- Optional tracking and logging of all activities performed on each sample
- Reports available as pdf or CSV files



Sample tracking report

### Tecan – Who we are

Tecan ([www.tecan.com](http://www.tecan.com)) is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics and clinical diagnostics. The company specializes in the development, production and distribution of automated workflow solutions for laboratories in the life sciences sector. Its clients include pharmaceutical and biotechnology companies, university research departments, forensic and diagnostic laboratories. As an original equipment manufacturer (OEM), Tecan is also a leader in developing and manufacturing OEM instruments and components that are then distributed by partner companies. Founded in Switzerland in 1980, the company has manufacturing, research and development sites in both Europe and North America and maintains a sales and service network in 52 countries.

**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **Denmark** +45 70 23 44 50 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170  
**Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Singapore** +65 644 41 886 **Spain** +34 935 95 2531 **Sweden** +46 31 75 44 000  
**Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33

Tecan Group Ltd. makes every effort to include accurate and up-to-date information, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided. Changes can be made at any time without notice. All mentioned trademarks are protected by law. For technical details and detailed procedures of the specifications provided please contact your Tecan representative. This may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

NucleoMag is a registered trademark of MACHERY-NAGEL GmbH & Co. KG, Germany. The NucleoMag kits developed designed distributed and sold by Macherey Nagel are intended for research use only.  
© 2014, Tecan Trading AG, Switzerland, all rights reserved.

[www.tecan.com/smartautomation](http://www.tecan.com/smartautomation)