



Automated nucleic acid purification using MACHERY-NAGEL NucleoSpin® kits

Freedom EVO® 75 optimized configuration

Automation doesn't get easier than this!

If you have limited lab space and wish hand-free, fully automated vacuum based isolation 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 food, feed, tissue, bacteria and cells. The solution offers flexible processing of batches single and multiple 8-well strips for up to 48 samples or 96-well plate for up to 96 samples.

Easy to use: Nucleic Acid Purification dedicated TouchTools™ 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



NucleoSpin principle



Nucleic acid purification workstation

Fast and efficient purification

Simplify your daily routine!

What can I use it for?

- Isolation of **genomic DNA from food and feed** with the NucleoSpin® 8 / 96 Food kit
- Isolation of **genomic DNA from human or animal tissue, cultured cells, or bacteria** with the NucleoSpin® 8 / 96 Tissue kit

What is included?

- Eight liquid handling channels
- Vacuum manifold
- Support for plates and disposable tips
- Robotic manipulator for automatic handling of the samples over the vacuum manifold
- 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.



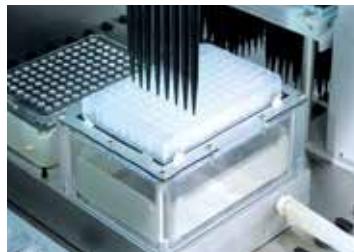
Nucleic acid purification wizard

Reliable optimized method

Use the 96-well binding plate for runs of 96 samples or 8-well strips for up to 48 samples.

Fast:

- 50 min for 48 samples
- 64 min for 96 samples



DNA purification by vacuum

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) 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.

NucleoSpin is a registered trademark of MACHERY-NAGEL GmbH & Co. KG, Germany. The NucleoSpin 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