
Service is a Tecan attribute

Service is a core competency of Tecan, ensuring that our customers get the most out of their investment. We strive to provide first-class care and support to our customers across the globe.

All services are provided by our broad network of experienced and certified field service specialists, guaranteeing the high level of service Tecan customers have come to expect.

Our services can be tailored according to your needs, and are provided through a range of Te-Care™ Service Contracts options or on demand.



Be Tecan trained and certified

We offer fast, convenient, effective training and certification for everyone who uses Tecan products.

Our training programs enable customers to build up the knowledge necessary to optimize the productivity of their Tecan instruments.

Our learning center provides online and offline training and is open to all, whether you want to learn the basics or gain more advanced tuition at the highest levels. Whatever your level of experience, we have a course that is right for you.



All data on page 5 was kindly provided by MACHEREY-NAGEL GmbH & CO. KG Neumann-Neander-Strasse 6; 52355 Düren, Germany.

Tecan are pioneers in automated liquid handling and innovative life science solutions. For over 30 years we continue to enable and support our customers to make the world a healthier and safer place.

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 48 174 **Portugal** +351 21 000 82 16 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74
Sweden +46 31 75 44 000 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **ROW** +41 44 922 8125

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this video or animation; 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 in this publication. Changes in this publication 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 in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at www.tecan.com/trademarks. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners.

© 2012, Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com



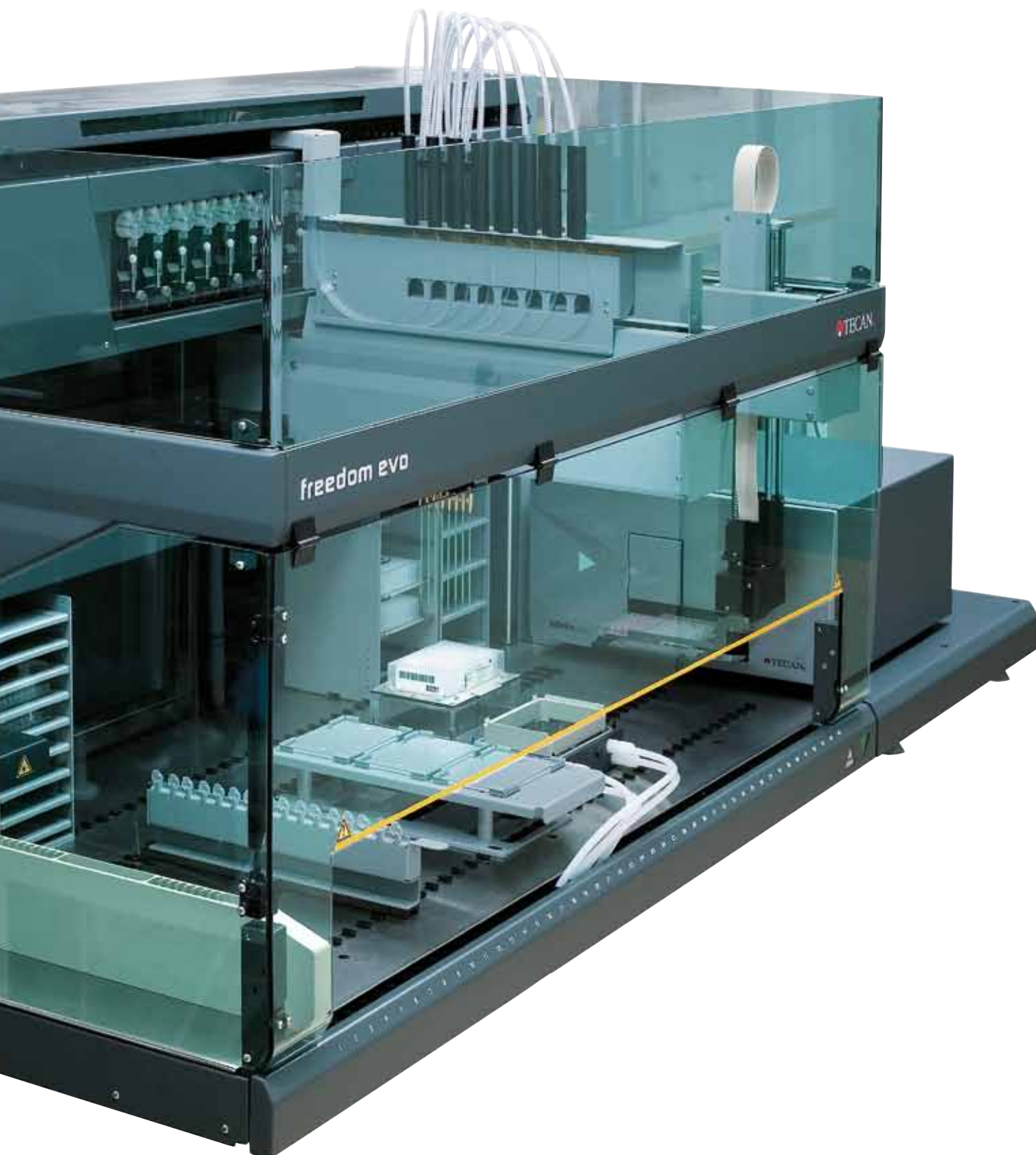


Accelerated nucleic acid purification

Flexible and proven solutions for Freedom EVO® nucleic acid purification

Need to free up time and increase productivity?

Tecan's Freedom EVO workstations are fully scalable, flexible systems that are ideally suited to nucleic acid purification and reaction set-up for a wide range of downstream applications.



Free choice of extraction chemistry

In collaboration with partners such as Promega and MACHEREY-NAGEL, we have established an extensive portfolio of co-developed and tested protocols for nucleic acid purification.



Choice of extraction methodology

Do you prefer solid phase extraction or magnetic bead separation? Our flexible Freedom EVO platforms allow both techniques to be performed on the same platform.



Flexible and upgradeable

A modular concept gives you the freedom to adjust your applications, meeting the changing demands of your workload and making the most of your laboratory's footprint and resources.



Fully integrated workflows

Purification, quantification and normalization of nucleic acids can be performed on the same platform, offering automation of your entire sample preparation workflow.



The Freedom EVO will address your nucleic acid purification needs of today and the future.

Strong partnerships to support your decisions

Tecan works closely with leading nucleic acid purification kit providers – such as Promega, Life Technologies and MACHEREY-NAGEL – to co-develop protocols for PCR and sequencing clean-up, as well as purification of genomic DNA, RNA, viral nucleic acids and plasmids from a wide range of samples, including:

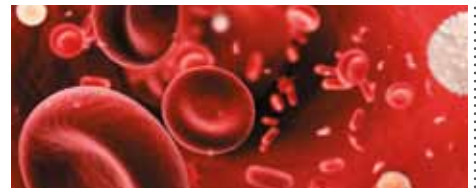
Plants

We meet your needs for fast and homogeneous extraction of high quality DNA from a variety of plants and fungi, from just a few milligrams up to 1.5 g of sample.



Blood

Both fresh and frozen human blood – and even more challenging samples such as animal blood – are purified with ease over an extensive volume range from 5 µl up to 10 ml.



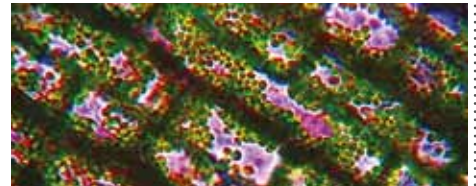
Body fluids

Viral nucleic acids can be purified in a true walkaway manner from up to 200 µl of body fluids, such as serum and plasma, for your large-scale screening projects.



Tissues

Flexible automated solutions for the isolation of genomic DNA from up to 20 mg of tissue material, without compromising on yield or purity.



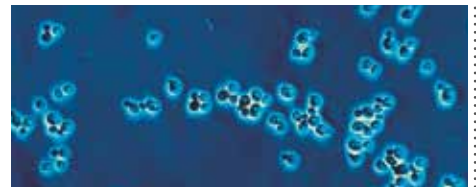
Cell cultures

Our automated platforms enable rapid, efficient and high quality purification of nucleic acid from up to 10^7 cells per sample.



Bacterial cultures

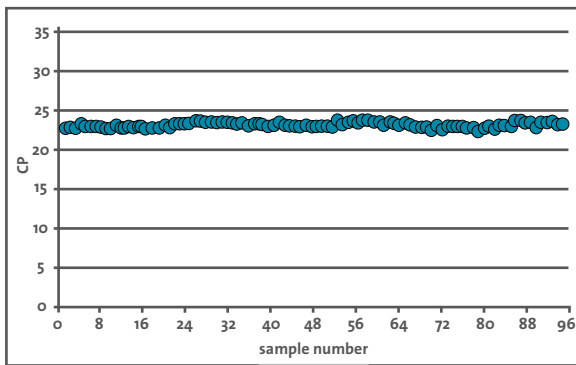
Pure genomic DNA or plasmids from up to 5 ml of your bacterial cultures, at high yields and speed.



Food products

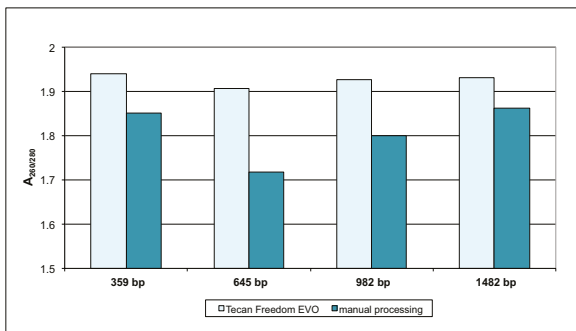
Pure nucleic acids from up to 200 mg of food or feed samples of plant or animal origin for your high throughput analyses.





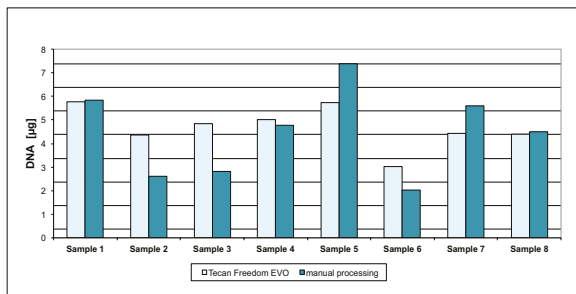
Highly reproducible results

Viral DNA extraction was performed with 96 identical samples and T7 DNA was determined using a real-time PCR system. Crossing points were highly reproducible at 23.3 ± 0.33 , with a CV of 1.42 %.



Highly pure nucleic acids

Each bar represents the average $A_{260/280}$ ratio from 24 extractions of plasmid DNA from 1.5 ml overnight bacteria cultures, processed either manually or using an automated method.



High yields

DNA yields from individual 200 µl samples of human whole blood range from 2 to 7 µg.

Choose your extraction chemistry

Solid phase extraction

For high throughput laboratories, we recommend our Te-VacS™ vacuum separation module. Each Te-VacS can accommodate 96-well filter plates, and the vacuum can be gradually increased to achieve optimal binding and elution conditions. Multiple Te-VacS units can be integrated onto one Freedom EVO platform to reach the desired sample throughput.

Magnetic bead extraction

For customers who prefer to work with a magnetic bead-based kit, a gripper arm is used to move plates to and from a static magnet, enabling high throughput applications and allowing almost any kit to be automated. For lower throughput applications, we offer our magnetic separation module, the Te-MagS™, which actively washes magnetic beads.

To view more data and download the latest application notes, please visit www.tecan.com/nap

The optimal solution for your needs

The unparalleled modularity of Freedom EVO workstations allow your liquid handling platform to be designed to meet your exact needs. Our specialists will be happy to discuss your optimal solution in full detail.

Enhance the capabilities of your Freedom EVO workstation

Third-party device integration

To extend your workflow beyond extraction, we can integrate almost any third-party device, such as thermocyclers and sealers, to complete your PCR set-up.

Quantification and normalization

Your complete nucleic acid workflow can be performed by installing a Tecan Infinite® series microplate reader beside or underneath the worktable, seamlessly integrating quantification and subsequent normalization after the extraction step.

Sample tracking and logistics

Automated logistics devices, such as hotels, a Carousel™ and Te-Stack™ modules, offer efficient and affordable ways to store and retrieve your samples. Process security is provided by integrated barcode reader options, or by the inclusion of a positive identification system (PosID™).

Shaking

To enhance the efficiency of your nucleic acid extraction protocol, the Te-Shake™ orbital shaker can vigorously shake most types of labware, including microplates, deep-well plates, PCR plates and tubes. It can also be configured or upgraded with different heating options to increase the yield during elution steps.

Centrifugation

Centrifugation steps are often crucial for sample preparation and the separation of individual layers prior to nucleic acid purification. The Freedom EVO portfolio accommodates a centrifuge beneath the worktable, maximizing the available deck space and minimizing the instrument footprint.



Sample tracking – PosID



Vacuum separation – TeVacS



Magnetic extraction – TeMagS



Magnetic extraction – 96-well magnetic separator



Shaking and heating



Thermocycling



Centrifugation



Quantification

Accessible automation

Easy-to-use software

Freedom EVOware's intuitive drag-and-drop graphical user interface (GUI) allows easy implementation of your methods. Process security is enhanced by Freedom EVOware's Safe Pathways feature, which reduces the risk of cross-contamination by guiding the tips around samples and other pieces of labware.



Seamless normalization

Freedom EVOware's Normalization Wizard enables all samples to be normalized to the same concentration. Dilutions can be calculated from data generated on integrated readers or by importing data files.



Streamlined sample tracking

Freedom EVOware's process security features enable sample tracking throughout the entire automated process. The built-in Sample Tracking module tracks and reports on samples for simplified LIMS integration, while password protection, audit trails and checksum help to detect and prevent unauthorized access.



Fast access

An optional touchscreen and TouchTools software increase the speed of access to previously written methods, further simplifying operation of the Freedom EVO platform.



Scheduling at your fingertips

The optional dynamic scheduler optimizes throughput for varying workloads, ensuring consistent processing, and increasing walkaway time.

