



Freedom EVO[®]

NGS workstation.

NGS LIBRARY PREPARATION & PCR SET-UP AT THE TOUCH OF A BUTTON

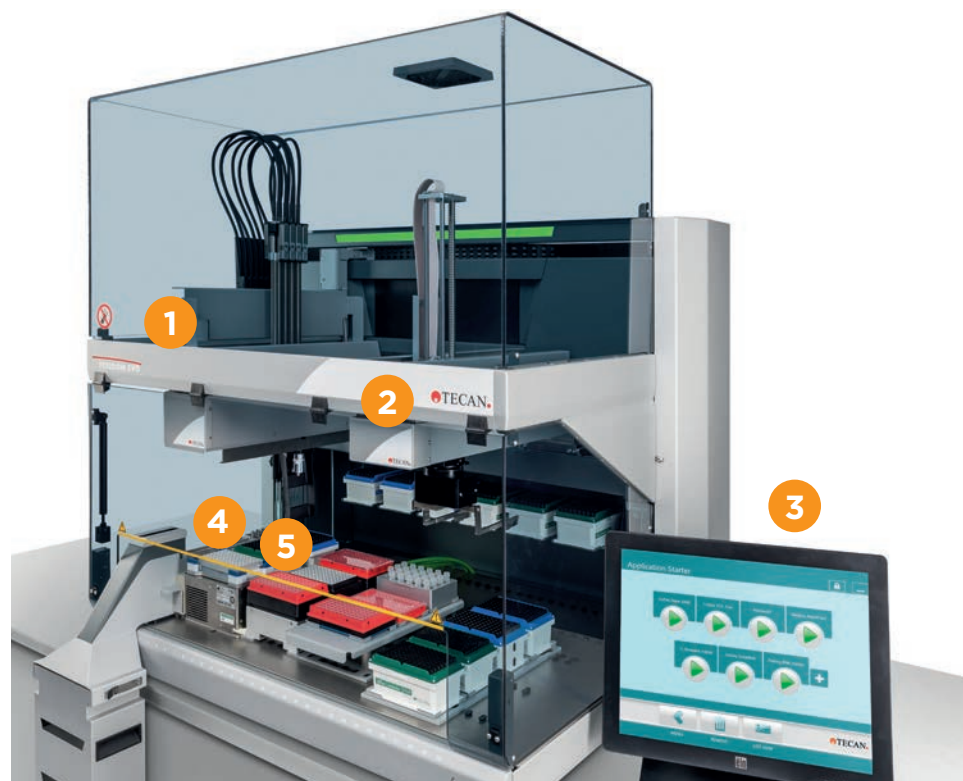


Next generation sequencing technology.

Next generation sequencing (NGS) technology continues to advance at an unprecedented speed and a growing range of new applications are being developed.

In conjunction with the ever increasing throughput of NGS instruments and the option to multiplex more and more samples, sample preparation has become a bottleneck. Automation of NGS sample preparation helps to reduce the workload, process more samples in parallel and increase the reproducibility of library preparations, while minimizing sample processing errors.

The Freedom EVO NGS workstation



Qualified protocols

Tecan developed and optimized automated NGS sample preparation protocols on the Freedom EVO NGS workstation in collaboration with key research centers and partners including Illumina® and IonTorrent™ to qualify all released protocols with sequencing data.

- 1 8-channel pipetting arm, (AirLiHa)
- 2 Robotic Manipulator Arm (RoMa)
- 3 Freedom EVOware® and TouchTools Suite™ software
- 4 Temperature controlled shaker
- 5 3 temperature controlled devices

Additional items include:

- Carrier for troughs, tip boxes, and PCR plates
- 96S Super Magnet
- Dust cover

The Freedom EVO NGS workstation, is based on the compact Freedom EVO 100 platform, equipped with the innovative air displacement pipetting technology. 8 pipetting channels allow highly precise pipetting from 1000 μ l down to 0.5 μ l. Variable tip spacing from 9 to 38 mm ensures efficient sample and reagent transfer from tubes to plates. The standard worktable layout also includes 3 temperature controlled units for reagent cooling and incubation of enzymatic reactions, as well as a shaker for efficient mixing of beads ensuring uniform yields. In addition, a simple touchscreen user interface makes protocol selection, instrument set-up and run fast and efficient.

What NGS sample preparation protocols are available?

Standard configuration workflows:

Illumina sample preparation:

- TruSeq® DNA Nano
- Truseq DNA PCR Free
- Nextera® XT
- Truseq Stranded mRNA
- Truseq Stranded Total RNA
- TruSeq RNA Access
- TruSight HLA v 2.0
- NEBNext Ultra II DNA

Capture protocols:

- Illumina Nextera Rapid Capture
- Agilent SureSelect XT™
- Roche SeqCap™ EZ

Ion Torrent sequencing:

- Ion AmpliSeq™
- Ion AmpliSeq panels



How long does it take to train an operator to use the instrument?

With the TouchTools™ graphic user interface, log in, selection of a protocol, worktable set up and starting a run requires minimal training and the user is guided through each step of the workflow.

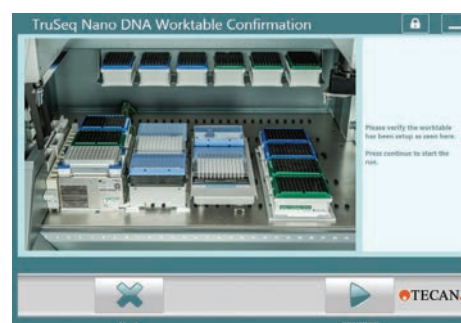


How is a protocol started?

The TouchTools application starter screen lets the operator select the desired protocol with the touch of a button.

Following protocol selection, specific run information including any variable number of samples from 1 to 96, final elution volume and other user defined parameters are entered.

Step by step user instructions avoid potential errors during worktable set-up to ensure correct initial conditions. Then a selection confirmation screen provides a run summary as a final check before the start of the run.



How does Tecan help customers to get started?

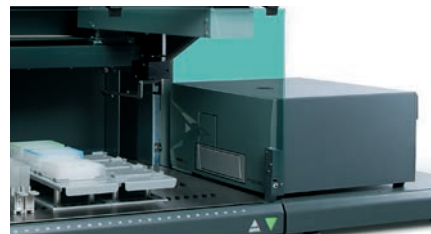
The Freedom EVO NGS workstation includes all currently available standard protocols at the time of installation. The installation process will include testing and training of one of the available protocols as agreed with the customer.

Custom script writing services are available for additional protocols, please ask your local Tecan sales representative for more information about this service.

ADDITIONAL OPTIONS:

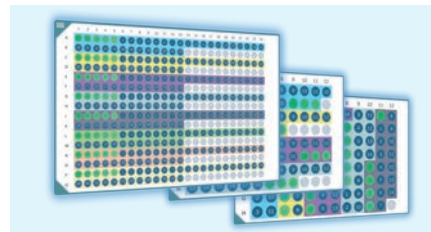
Fully automated quantification and normalization

- Integration of Infinite F200 reader & Normalization Wizard



TouchTools™ PCR Wizard

- Effortless PCR set-up for your everyday needs



CONSUMABLES:

- 30057813 Disposable Tips, 50 µl, Conductive, Filtered, Tecan Pure Purity Level, 24 x 96 pcs, for refill in disposable tips SBS box
- 30057815 Disposable Tips, 200 µl, Conductive, Filtered, Tecan Pure Purity Level, 24 x 96 pcs, for refill in disposable tips SBS box
- 30058506 Disposable Tips refill box, SBS format, 10 pcs
- 30055743 Disposable Troughs, 25 mL, PP, grey, Tecan Pure Purity Level, 120 pcs
- 10613049 Disposable Troughs, 100 mL, PP, grey, Tecan Standard Purity Level, 108 pcs

SERVICE AND TRAINING ARE A TECAN ATTRIBUTE

Tecan's service and training products are designed to ensure your systems's productivity. Our global network of certified field service specialists delivers a high level of support when you need it. Services are provided through a range of TeCare™ Service Contract Options or on demand.

Empowering users through training ensures maximum productivity from Tecan products. Tecan delivers trainings to meet the needs of diverse users through classroom, onsite and online settings. For more information, please visit www.tecan.com or contact your local Tecan sales representative.

.....
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 93 490 01 74
Sweden +46 8 750 39 40 **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 within this publication, 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. 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 <http://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. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland.
© 2016 Tecan Trading AG, Switzerland, all rights reserved.

For research use only. Not for use in diagnostic procedures.

www.tecan.com



