

VEYA

EFFORTLESS AUTOMATION.



EFFORTLESS AUTOMATION.

Veya™ is a multiomics platform designed to automate both research and clinical workflows. It combines digital tools to streamline operations, offering an intuitive experience that makes automation easy. With predeveloped scripts for quick and easy setup, critical tasks can be automated effortlessly. Backed by Tecan's proven technology, Veya delivers reliable performance across various lab environments.



Effortless set-up, with predeveloped protocols¹ and an intuitive user interface.



Fewer operating errors thanks to cutting-edge digital tools and automated environmental calibration.



Optimized productivity and up to 64 % higher instrument utilization with intelligent analytics and OneView™ display.



Personalized, digitally empowered service at every step of the customer journey, with a 24-hour remote response time².



Excellence in regulated environments, including CE-IVD³ compliance.



Powered by Tecan's trusted modular platform, proven in clinical markets, with validated performance when used with Tecan consumables.



¹ Protocols need to be validated by the user.

² For advanced and premium service contracts.

³ Compliant to IVDR (Regulation (EU) 2017/746) class A; USA Class 1 IVD, 510(k) exempt.

EFFORTLESS MULTIOMICS.

Veya multiomic suites are specialized configurations with ready-to-use predeveloped scripts⁵, designed to streamline the automation of complex multiomic workflows. Whether scaling up or adjusting workflows to meet specific needs, vControl™ simplifies method creation with its intuitive interface and Smart Commands.

Veya NAP Suite

- Automation for a wide variety of magnetic bead-based nucleic acid purification protocols.
- Walk-away solution for up to 96 samples in less than three hours.*
- Standardized worktable and wide volume range pipetting to track and extract samples at ease.
- Enhanced genomic workflows with safe sample transfer and PCR setup.

Veya NGS Suite

- Walk-away solution for any number of samples, up to 48, processed in just four hours – without the limitations of fixed sample increments.*
- Universal worktable, compatible with short- and long-read protocols and any type of DNA or RNA sample.
- Future-proof with an extensive and growing applications library, compatible with Fluent® Automation Workstation.

Veya Cell Biology Suite

- Sterile cell culture environment, with the integrated HEPA filter⁶.
- Streamlined cell-based assays using built-in smart commands for efficiency and reliability.

Veya Clinical Suite

- Ideal entry-level solution for clinical workflows, designed for sample distribution to various daughter plates.

**Available methods
for NAP and NGS**



* The required time is determined by the protocol.

⁵ Protocols need to be validated by the user.

⁶ The cell-based experiments were conducted under RUO conditions.

EFFORTLESS OPERATION.

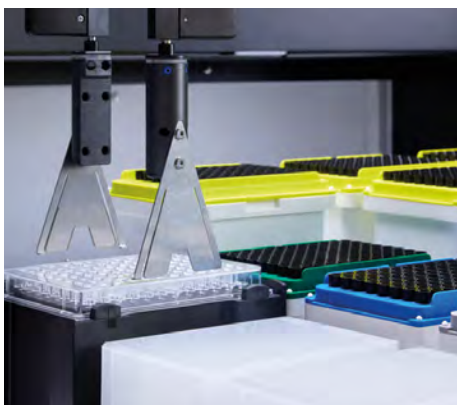
Veya features a unique auto-adjusting liquid handling technology with a versatile 1 μ l to 5 ml pipetting range and customizable gripper fingers. With up to 34 plate positions and additional modules for cooling, heating, shaking and decontamination, this system is powered by robust software and AI tools for performance, quality and ease of use.





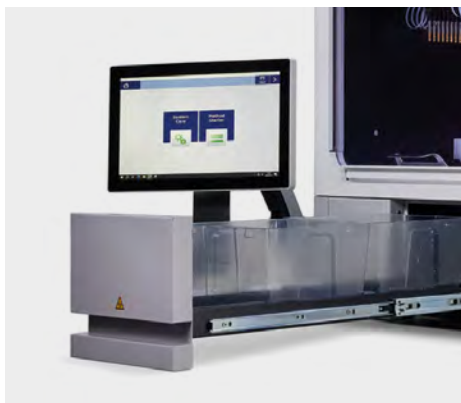
Plate handling.

Easy, on-deck labware handling with the Channel Gripper. Access areas below the deck with an optional robotic gripper arm, without consuming space on deck.



Tecan tips only.

A wide range of disposable tips are available for use with Veya. Only original Tecan tips, designed and validated alongside our instruments, ensure reliable and reproducible performance.



Waste management.

The tip-eject system prevents cross-contamination by using a closed ejection station and a sealed waste drawer beneath the deck. Waste bins can be autoclaved for easy sterilization.



Touch screen.

A convenient way for operators to interact with vControl software.



EFFORTLESS USER EXPERIENCE.

Veya provides real-time insights into system status and guides operators through every step. It automatically detects loading and pipetting errors, while advanced analytics monitor instrument performance and can increase utilization by up to 64%.



OneView.

Provides real-time data on system status:

- Instrument system care and environmental status.
- Real-time run status and history of last runs.
- Introspect™ efficiency score.
- Direct access to Tecan's self-help knowledge portal, available 24 hours a day.



Introspect.

Tecan's advanced data analytics software provides actionable insights to increase instrument utilization and performance, along with remote monitoring capabilities for extended walkaway times.



LabNavigator™.

Intuitive software guides lab operators through end-to-end workflows using a single, intuitive interface (optional).



DeckCheck™.

An integral sensor system that detects loading errors, ensuring smooth operations and successful runs.

Pressure Monitored Pipetting.

Continuous AI-based monitoring of pipetting pressure identifies pipetting errors in real time.

EFFORTLESS SUPPORT.

Every Veya comes with access to standard support services, including self-help tools and basic e-learning resources. These services are designed to give users quick, easy access to troubleshooting and ensure smooth operations.

Personalized service and digital support.

For more comprehensive support, a range of Te-Care™ service contracts is available. All service contracts include remote instrument diagnostics and support services, ensuring timely responses, with a remote response commitment within 24 hours⁷, minimizing downtime and keeping Veya at peak performance.

Explore the available service contracts to ensure that your lab receives the care and support that it needs.

Overview of service contracts.



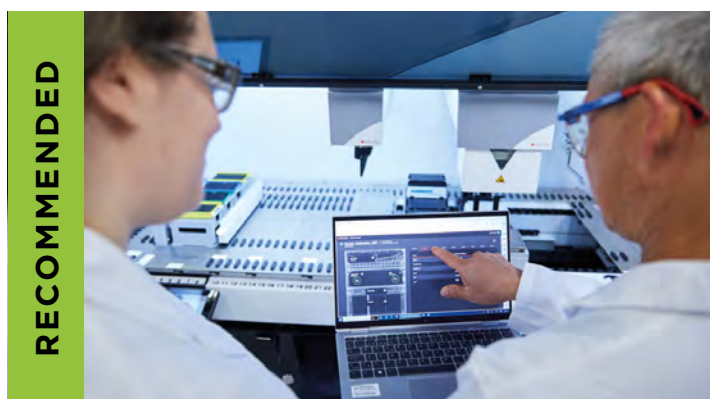
Te-Care CORE.

Efficient remote troubleshooting with timely access to Tecan specialists for quick support.



Te-Care BASIC.

Optimal performance through regular onsite preventive maintenance conducted by qualified and certified service personnel.



Te-Care ADVANCED.

High performance with onsite preventive maintenance, remote troubleshooting before repairs and prioritized support.

Te-Care PREMIUM.

Peace of mind with onsite preventive maintenance, complete onsite troubleshooting and priority support.

⁷ For Te-Care ADVANCED and PREMIUM service contracts.

TECHNICAL FEATURES AND SPECIFICATIONS.

General

1 OneView

2 Status light

3 4 or 8 channels

4 1 µl – 5 ml

5 Air Restriction Pipetting (ARP)

6 DeckCheck

+ Also available as an IVD medical device.

7 vControl

8 Modular worktable

9 Two sizes: up to 22 to 34 plate positions/6 and 9 segments

10 Introspect

Ready-to-use application suites

Cell Biology	NAP	NGS	Clinical*
--------------	-----	-----	-----------

* reagent kits available on request

For further options, please contact your local Tecan sales representative.

LEARN MORE

Contact Tecan sales organization for local regulatory status of the product.

Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **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 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **South Korea** +82 2 818 3301
Spain +34 93 595 25 31 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 408 782 5600 **Other countries** +41 44 922 81 11

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.

© 2025 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com

TECAN

403123 V1.0, 2025-01