

Resolvex® A200 Omics and Resolvex A200 24.

THE NEXT GENERATION OF RESOLVEX A200 POSITIVE PRESSURE PROCESSORS



Resolvex A200.

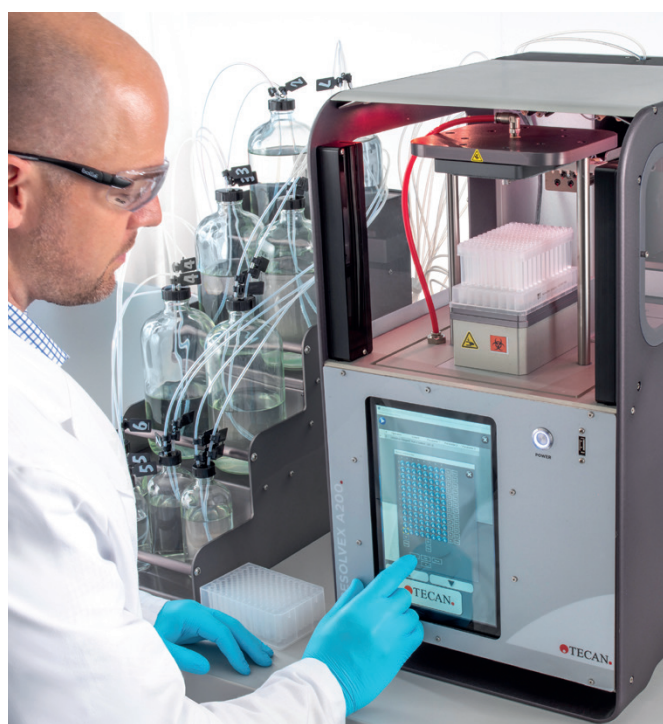
BENCHTOP, AUTOMATED POSITIVE PRESSURE PROCESSOR

The Resolvex A200 is an industry-proven, standalone system for automated positive pressure sample preparation. This compact processing solution provides precise dispensing of up to 11 solvents – from water to organic solvents – using industry-leading programmable pressure profiles to optimize sample loading and improve accuracy, precision and method robustness.

The next generation Resolvex A200 Omics and Resolvex A200 24 builds on this success, providing additional features that enable scientists to significantly expand the platform's workflow capabilities.

The Resolvex A200 Omics incorporates a HEPA filter air system to reduce particulates, enabling applications sensitive to air contamination. Particulates – for example, keratin – can impact the quality of results when working with mass spectrometry (MS). This is of particular concern with biochemical and molecular biology MS workflows.

The Resolvex A200 24 gives users the capability to use larger format sample prep consumables, including 3 and 6 ml columns, and 24-well plates. Cleverly designed accessories allow users to run additional workflows, such as large-scale, giga-sized nucleic acid purification (NAP) methods, improving throughput and enhancing method robustness by reducing the potential for human errors.



Applications:

- Bioanalysis
- Analyte clean-up and enrichment for therapeutic drug monitoring, drug screening and other clinical chemistry workflows
- Nucleic acid screening and purification
- Life sciences research
- Food safety testing
- Environmental monitoring

ResolveX A200 Omics

Features and benefits

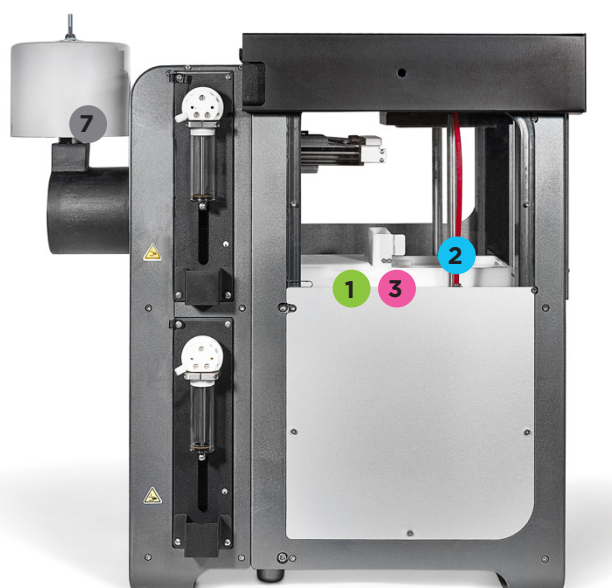
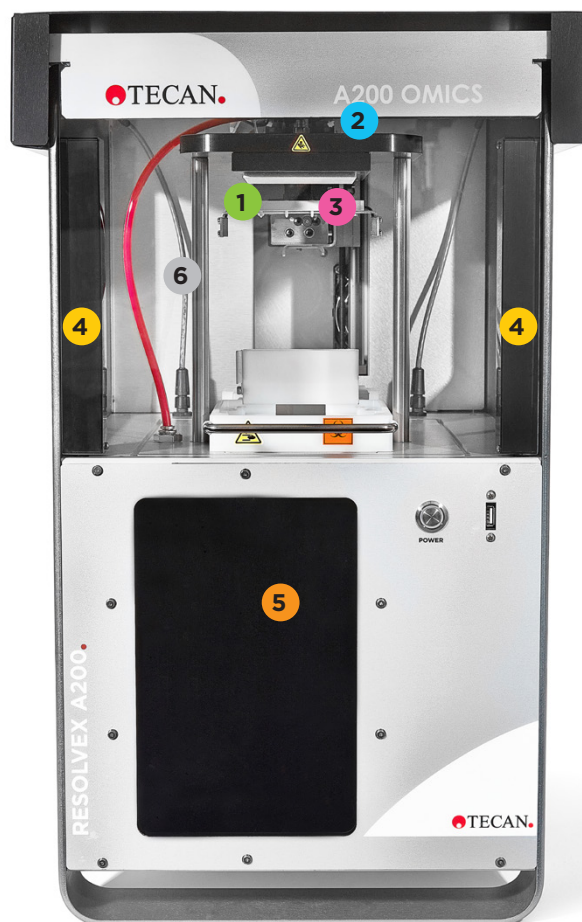
1. Sensor-controlled movement to perfectly position the dispense head at each step

3. Automated liquid dispenser

- Tecan Cavo® dual syringe pump increases dispense speeds for high volume sample preparation processes
- Accurate delivery of protocol solvents
- Connection of up to 11 solvents at a time
- Motion-controlled 8- or 4-port dispense head

2. 24- or 96-channel positive pressure manifold

- Use of positive gas pressure provides maximum reproducibility
- Uniform pressure processing across columns and plate wells
- Pneumatic clamping system with a dual air flow path ensures consistent and reproducible pressures
- Precise flow control reduces processing times and enhances performance



4. Safety light curtain detects presence of objects to prevent possible damage or injury

6. Enclosed magnetic cover to prevent particles from entering the unit

7. HEPA filter delivers consistent filtered air into the unit, significantly reducing particulates

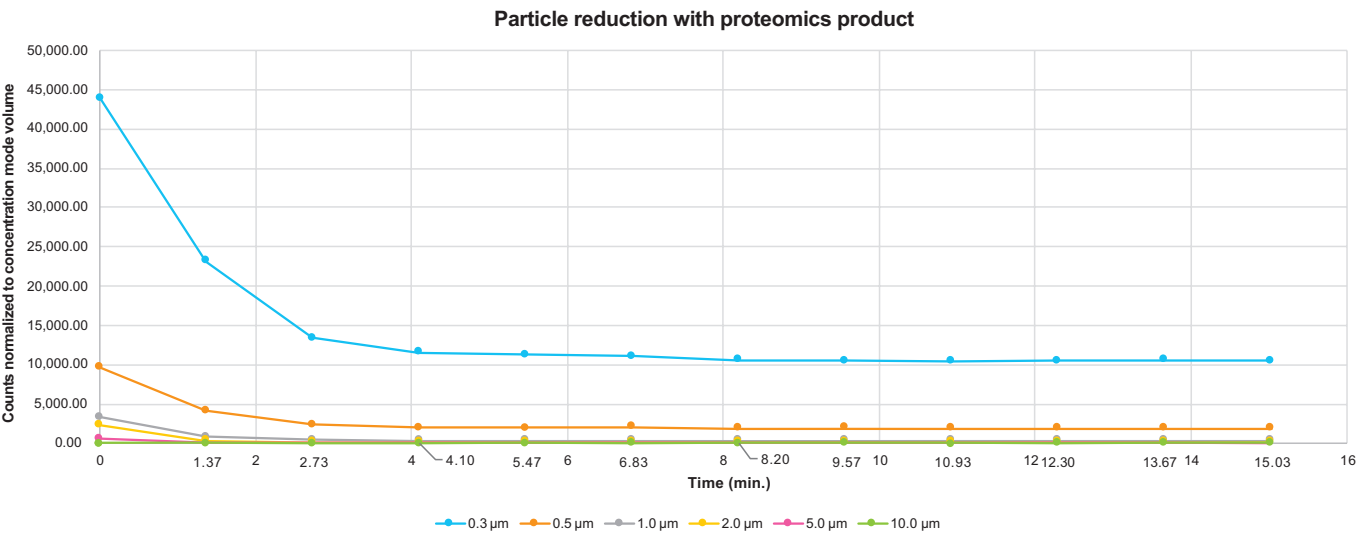
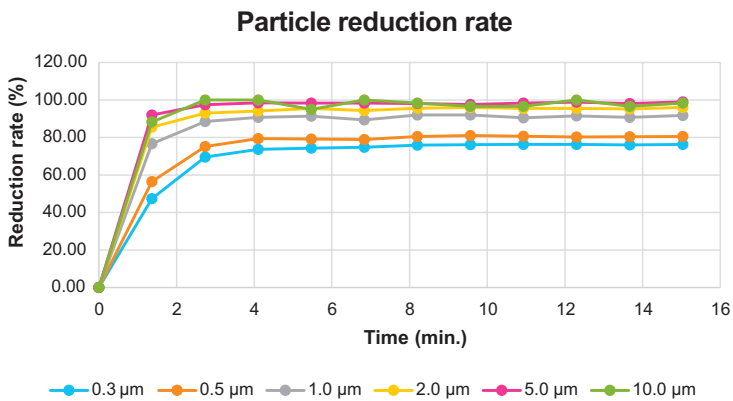
5.

- Microsoft Windows® PC application and intuitive touchscreen interface
- Programmable solvent dispensing
- Precise pressure profiles for accurate and optimized flow control
- Ability to program and run multiple methods on a single system



Resolvex A200 Omics HEPA performance

The Resolvex A200 Omics provides an enclosed environment to significantly reduce airborne particulates and ensure a constant airflow in the processor workspace, minimizing the possibility of sample contamination. As a standalone unit, the Resolvex A200 Omics can be operated anywhere in the lab, while still providing the correct sample preparation environment.



ResolveX A200 24

Features and benefits

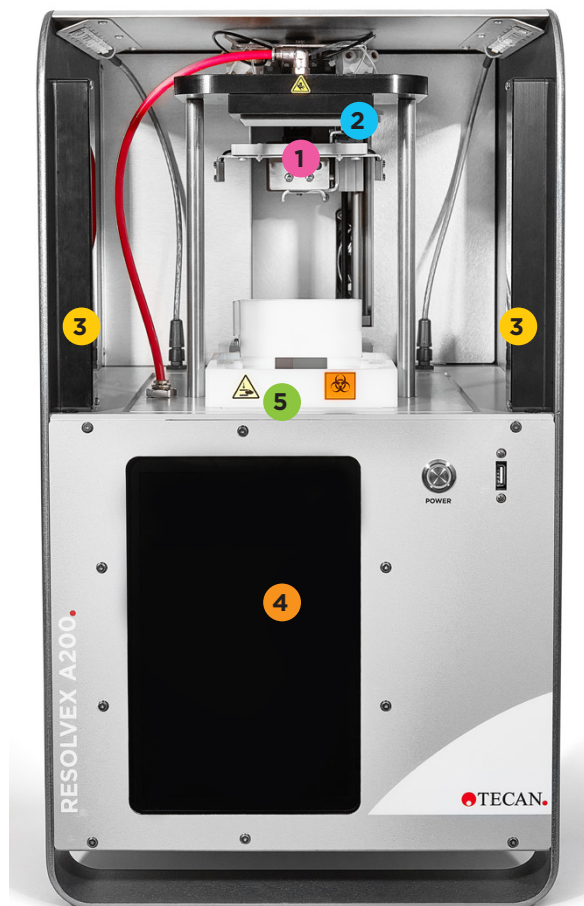
1. Automated liquid dispenser

- Tecan Cattro dual syringe pump increases dispense speeds for high volume sample preparation processes
- Accurate delivery of protocol solvents
- Connection of up to 11 solvents at a time
- Motion-controlled 4-port dispense head

- ### 3. Safety light curtain
- detects presence of objects to prevent possible damage or injury

4.

- Microsoft Windows PC application and intuitive touchscreen interface
- Programmable solvent dispensing
- Precise pressure profiles for accurate and optimized flow control
- Ability to program and run multiple methods on a single system



2. 24-channel positive pressure manifold

- Use of positive gas pressure provides maximum reproducibility
- Uniform pressure processing across columns and plate wells
- Pneumatic clamping system with a dual air flow path ensures consistent and reproducible pressures
- Precise flow control reduces processing times and enhances performance

5. Updated and redesigned waste station

- Easier to use and replaceable to reduce long-term contamination
- Nest design to allow for easier placement of plates and racks/spacers
- Lower profile to accommodate larger columns and stacks
- Solid, high density plastic to reduce contamination
- More automation friendly for robotic transfer of samples

Resolvex A200 24.

ADDITIONAL FEATURES



Flexibility to process from one to 96 samples at the touch of a button



3 ml



6 ml

Single tube racks for 1, 3, and 6 ml columns offer flexibility and avoid waste

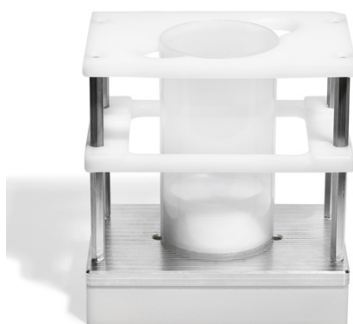


16 x 100 mm test tube and auto sample injection vial racks for different elution needs

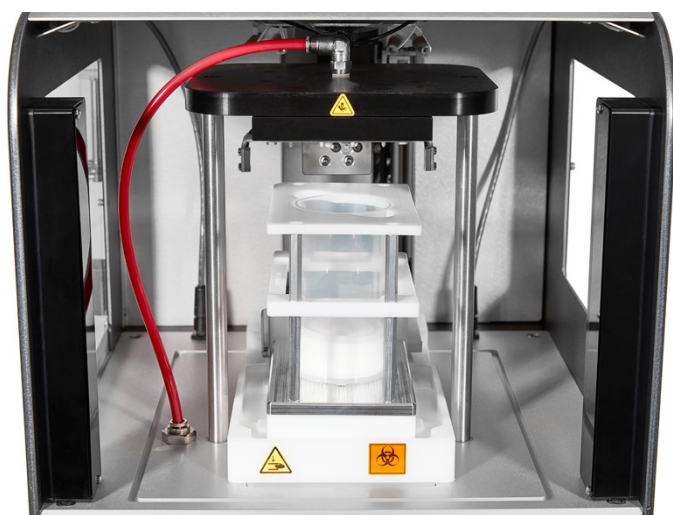


Resolvex A200 24.

ADDITIONAL WORKFLOW ADAPTERS



Specially designed giga column adapter racks for the MACHEREY-NAGEL Giga column system



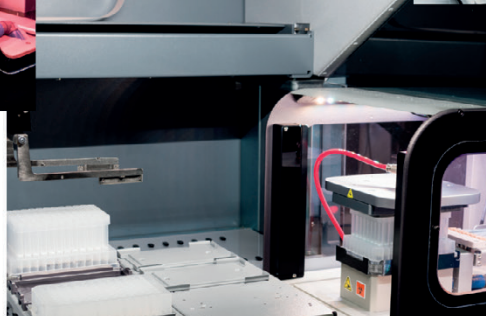
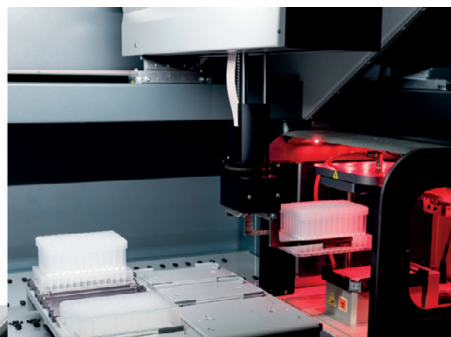
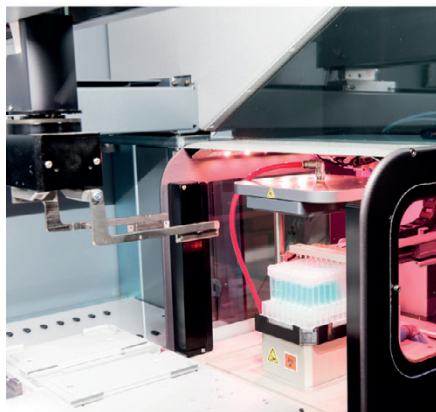
Specially designed focus adapter to prevent splashing and possible contamination

MACHEREY-NAGEL Giga column set-up on the Resolvex A200 24 system

A complete sample handling and positive pressure processing workflow solution

Resolvex A200 integrated with a Fluent® or Freedom EVO® liquid handling platform

Combining Tecan's proven liquid handling platforms with the industry-leading Resolvex A200 positive pressure processor provides a fully automated pre-analytical sample processing workstation.



+1 626 962 0010

customerservice-spe@tecan.com

www.tecan.com/resolvex-a200-next-generation-positive-pressure-processing

For Research Use Only – not for use in diagnostic procedures.

.....
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 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **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.
© 2023 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com

