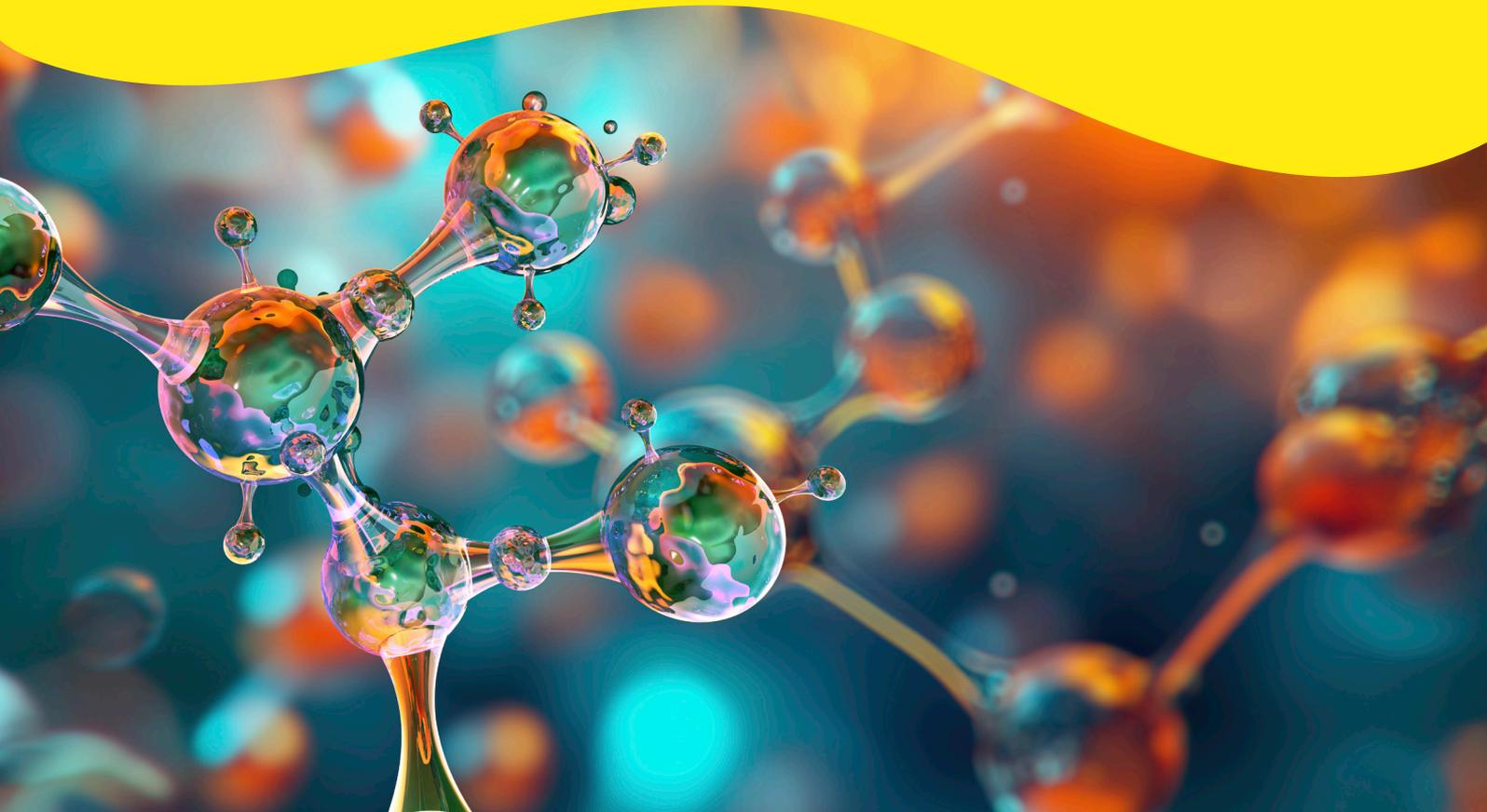


revvity

LIFE SCIENCE RESEARCH

Automated liquid handling solutions: from benchtop to integration.

Comprehensive guide



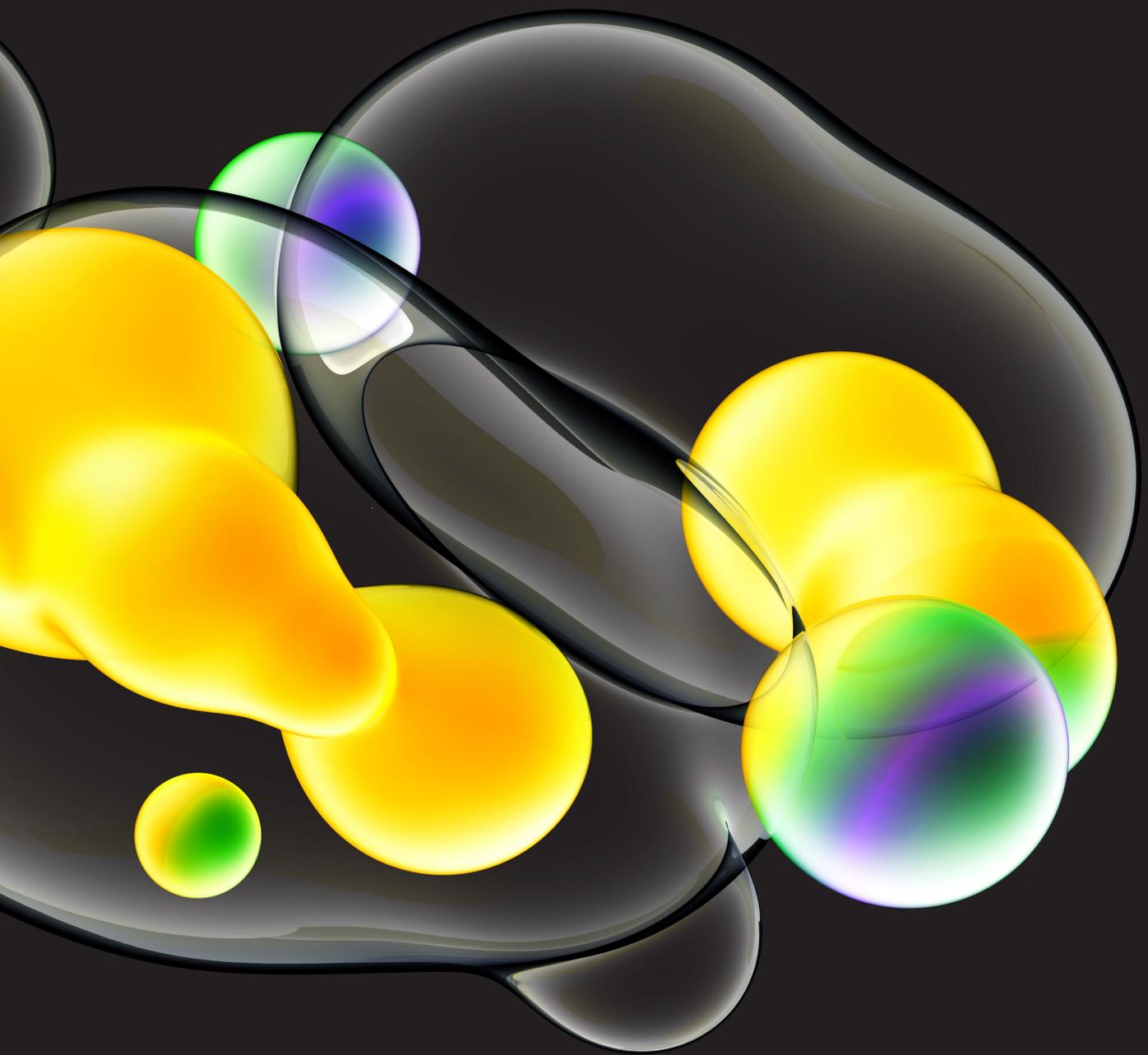


Table of contents

Introduction	3
Section 1 - Automated liquid handling portfolio	3
AssayMate™ workstation: The Intelligent Automation Entry Point.	4
Zephyr™ G3 workstation – Robust Mid-Range Workstation.	5
Fontus™ and JANUS™ workstations – The Flexible Powerhouses	6
The Fontus™ Workstation: Next-Generation Adaptability	6
The JANUS™ G3 Workstation: The Versatile Performer	8
Fontus and JANUS systems’ technical comparison	9
FlexDrop™ Plus dispenser – contactless nanoliter precision	10
Section 2: The Science of the Dispense	11
Revvity Pipetting Technology Decision Matrix	12
Section 3: Applications, the solution map	12
Guide by application	13
Guide by sample type	13
Application-Dedicated Workstations	14
Section 4: Better Together – platform synergy	15
Section 5: explorer™ G3 workstation - Integrated Automated Solutions	17
Section 6: Why Revvity – Your Partner in Discovery	18
Conclusion: Challenge Accepted	20

Introduction

In today’s fast-paced research environment, the bottleneck to breakthrough discovery is rarely the lack of ideas but rather the limitation of manual processes or difficult repeatability. Repetitive pipetting, variability in liquid handling, and the increasing pressure to miniaturize assays without sacrificing data quality could slow the momentum of even the most innovative labs without proper equipment.

We believe that automation should be accessible, scalable, and, above all, scientifically driven. Whether you are a startup looking for your first benchtop workstation or a larger laboratory requiring complex, 24/7 integrated workflows, our portfolio of solutions helps offer the harmony of precision, speed, and adaptability.

By helping you choose the liquid handler that fits your application needs, you will be able to support your goal to reduce human error and reagent waste, and achieve the reproducibility required for the next generation of therapeutics.

Revvity’s liquid handling portfolio helps to return time to the scientist.

Welcome to a more productive lab.

Welcome to Revvity.

SECTION 1

Automated liquid handling portfolio

Flexibility without sacrificing precision

Automating your science, from benchtop to fully automated workflow

Revvity's automated liquid handling solutions are built on the philosophy of providing workflow flexibility to minimize errors, reduce hands-on time, and increase reproducibility without overcomplicating the user experience.

Whether you are performing delicate sub-microliter dispensing or managing high-volume sample preparation, our portfolio offers solutions to automate your science at any stage of laboratory growth.

From the compact, affordable AssayMate™ workstation designed for any benchtop, to the high-throughput, larger scale explorer™ G3 workstation integrated solutions, Revvity systems offer adaptability in throughput, capacity, and dynamic volume range. Our workstations are more than just robots; they are comprehensive application-focused solutions equipped with easy-to-use software, verified protocols, and remote workflow monitoring to enable your lab to stay productive and future-ready.

Explore our portfolio below to find the system that matches your current throughput needs and can help you scale your future discoveries.



[Learn more](#)

Scan or click
to Learn more!

AssayMate workstation: The intelligent automation entry point

Compact and intelligent: The AssayMate™ liquid handler is a benchtop-sized handler that bridges the gap between manual pipetting and robotics. Features pressure-based liquid level detection (PLLD) and a reconfigurable recessed deck for heater/shaker modules.

- **Compact performance:** A benchtop-sized handler designed to help bridge the gap between manual pipetting and full-scale robotics.
- **Wide volume range:** Handles up to 1000 µL without needing to switch heads, supporting a fixed 8-channel head and gripper configuration.
- **Intelligent features:** Integrates pressure-based liquid level detection (PLLD) for reliable operation. Workflow error handling mechanism*, such as detecting a missing plate lid when the protocol calls for it, allows for quick experiment resumption.
- **Modular deck:** Features a reconfigurable recessed section for optional modules like heater shakers, heater coolers, and magnetic blocks.
- **User-friendly:** Managed by the AssayPREP™ software with a simple drag-and-drop interface requiring minimal training.



YAS8G01

* For most error conditions.

Resources:

- [AssayMate workstation brochure](#)
- [AssayMate workstation specification sheet](#)
- [AssayMate technical overview webinar \(incl. a short software demo\)](#)



Learn more

More information
and resources

Zephyr G3 workstation – robust mid-range workstation

The Zephyr™ G3 workstation is a compact, cost-effective, multi-channel liquid handler.

- **Simple and robust:** A compact, 12-position workstation designed to support reliability.
- **High-volume efficiency:** Equipped with a 96-channel high-volume head (HVH) capable of processing an entire plate in minutes with high precision (CV <2% for 5-200 µL).
- **Contamination control:** Features an innovative offset gripper that holds lids while pipetting, helping prevent deck-surface contamination.
- **Versatility:** Supports partial-tip loading (single rows/columns).
- **Extended capacity:** Seamless integration with the Twister™ III plate handler robotic arm increases storage capacity (tips, plates) and optional additional modules integration (e.g. ODTc).



Zephyr G3 Low-Volume Head (LVH) model (CLS145301) includes:

- 96/384-well LVH head
- Left gripper
- Tip array options: 96- or 384- mandrel array
- LVH head utilizes disposable tips
- Dynamic volume range: 1-25 µL

Zephyr G3 High-Volume Head (HVH) model (CLS145302) includes:

- 96-well HVH head
- Left gripper
- HVH head utilizes disposable tips
- Dynamic volume range: 1-200 µL

Resources:

- [Zephyr workstation specification sheet](#)
- [Twister III robotic arm specification sheet](#)
- [Bioz citations](#)



Learn more

More information
and resources

Fontus and JANUS workstations – the flexible powerhouses

The Fontus™ and JANUS™ platforms are Revvity's flagship workstations for labs that need the combination of high-capacity throughput and intricate, variable-volume pipetting.

These systems offer modular deck sizes, Mini, Standard, and Expanded, as well as an Integrator model for the JANUS workstation. They also offer two pipetting technologies with the 96- or 384-channel main array head and the Varispan™ 8-span head for advanced tubes to plates reformatting capabilities.

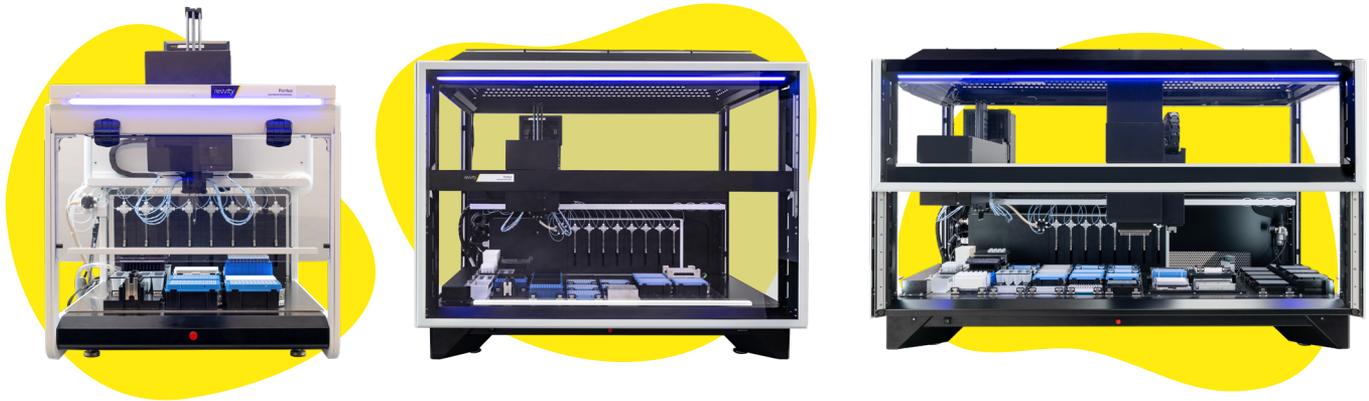
Fontus and JANUS workstations support FDA 21 CFR Part 11 compliance. The WinPREP™ software contains built-in technical controls and features specifically designed to support FDA 21 CFR Part 11 regulations.

The Fontus and JANUS workstations include both dedicated, application-specific workstations and personalized automated solutions.

The Fontus workstation: next-generation adaptability

The Fontus platform represents the evolution of Revvity's liquid handling, designed to “defy obsolescence” by scaling alongside your research.

- **Varispan™ pipetting head:** Equipped with 8 independent pipetting channels that move in the X, Y, and Z axes. This allows the transfer of samples between different labware types (e.g., from tubes to 384-well plates) in a single step.
- **Enhanced tip flexibility:** Features VersaTip™ technology, allowing the system to use a mix of fixed, washable tips and disposable tips within the same protocol—perfect for reducing cost-per-run without sacrificing speed.
- **Accusense™ technology:** Patented liquid level sensing that can detect even non-ionic polar solutions and solvents like DMSO to support reliable aspirations.
- **Intuitive user experience:** Powered by WinPREP™ software with an Application Assistant that simplifies method creation through a visual, drag-and-drop interface.
- **Enhanced ergonomics** with new enclosure and optional UV light with entire deck coverage
- **Modular growth path:** Start with a single-arm Mini or Standard deck and, as your throughput needs increase or your workflows become more complex, advance to a dual-arm Expanded configuration, transferring your protocols easily.
- **High-capacity storage:** Includes sunken tip box storage on the deck and the ability to link with PlateStak™ Automated Microplate Handler ([CLS156605](#)) for massive walkaway capacity.

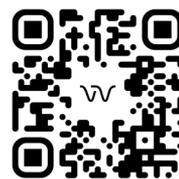


The Fontus workstation family includes the following models:

- Fontus 8i Mini Workstation, 8-Tip Varispan ([YFS8002](#))
- Fontus Workstation, Single-Arm, 8-Tip Varispan, Standard (YFM8001)
- Fontus Workstation, Single-Arm, 8-Tip Varispan, Expanded (YFL8001)
- Fontus Workstation, Dual-Arm, 8-Tip Varispan, 96 Main Array, Standard ([YFM8H01](#))
- Fontus Workstation, Dual-Arm, 8-Tip Varispan, 96 Main Array, Expanded ([YFL8H01](#))
- Fontus Workstation, Dual-Arm, 8-Tip Varispan, 384 Main Array, Standard ([YFM8L01](#))
- Fontus Workstation, Dual-Arm, 8-Tip Varispan, 384 Main Array, Expanded ([YFL8L01](#))

Resources:

- [Fontus 8i Mini workstation flyer](#)
- [Fontus workstations specification sheet](#)
- [Fontus workstation video](#)
- [PlateStak Plate Handler flyer](#)
- [Immunoassays workflow solutions infographic](#)
- [Fontus webinar on-demand: technical overview](#)



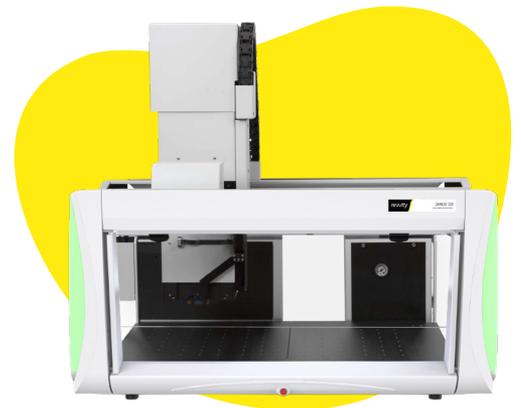
[Learn more](#)

**More information
and resources**

The JANUS G3 Workstation: The versatile performer

The JANUS™ G3 has become an industry standard for labs requiring an extensive dynamic volume range and precise movement across diverse labware.

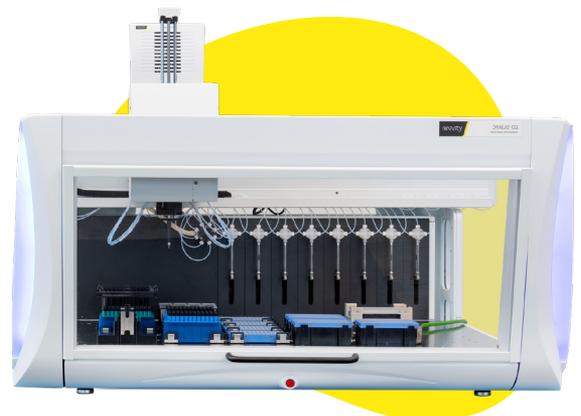
- **Varispan™ pipetting head:** Equipped with 8 independent pipetting channels that move in the X, Y, and Z axes. This allows the transfer of samples between different labware types (e.g., from tubes to 384-well plates) in a single step.
- **Accusense™ technology:** Patented liquid level sensing that can detect even non-ionic polar solutions and solvents like DMSO to support reliable aspirations.
- **Massive volume range:** Delivers consistent and reproducible results across a range from 0.5 µL to 5000 µL.
- **Additional capability:** Ability to link with PlateStak™ Automated Microplate Handler ([CLS156605](#)) for extended storage and walkaway capacity.
- **Application-specific specialization:** The JANUS architecture serves as the foundation for highly specialized workstations, such as the JANUS BioTx system for protein purification, which integrates vacuum manifolds and automated fraction collection.

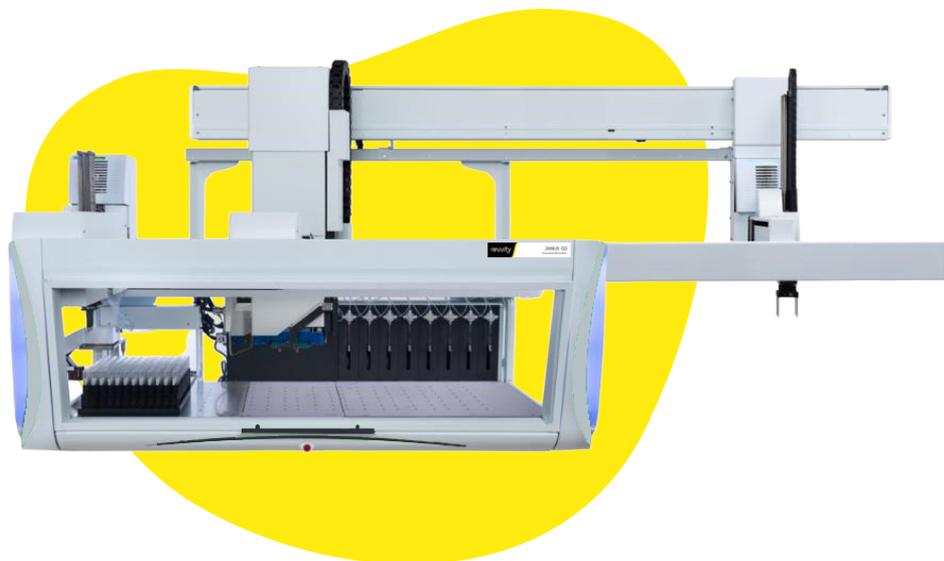


The JANUS workstation family includes the following models:

- JANUS G3 Mini Workstation, 8-Tip Varispan ([YJS8001](#))
- JANUS G3 MDT Workstation, single-arm MDT, expanded ([YJLM001](#))
- JANUS G3 MDT Workstation, single-arm MDT, expanded, with gripper ([YJLMG01](#))
- JANUS G3 Workstation, Dual-arm 8-Tip Varispan and MDT, standard ([YJM8M01](#))
- JANUS G3 Workstation, Dual-arm 8-Tip Varispan and MDT, expanded ([YJL8M01](#))

* Enclosure available for some models.





JANUS G3 Integrator model

Resources:

- [JANUS G3 workstations specification sheet](#)
- [AlphaLISA™ assay automation on JANUS workstation](#)
- [Bioz citations](#)



[Learn more](#)

More information
and resources

Fontus and JANUS systems' technical comparison

To help you navigate between Revvity's most versatile platforms, the following table summarizes the Fontus and JANUS G3 technical specifications and deck architectures.

Feature	Fontus workstation	JANUS G3 workstation
Configurations	2 arm configurations: single or dual arm	3 arm configurations: single, dual arm, or Integrator with additional arm for extra gripper
Available deck sizes	Mini, Standard and Expanded. (Mini available as Varispan-only)	Mini, Standard, Expanded, and Integrator.
Software interface	WinPREP™ with Application Assistant	WinPREP™ with Application Assistant
Pipetting technology	Liquid Displacement (Varispan) & Air Displacement (Main Array).	Liquid Displacement (Varispan) & Air Displacement (96/384 MDT) with software-controlled swap during workflow.
Mini deck capability	Varispan Only: Ideal for tube-to-plate transfers in a small footprint.	Varispan (span-8-channel) or MDT (96- or 384-channel)
Deck capacity	Higher: Standard: 28+4 / Expanded: 36+4. Optimized for high-density layouts.	Standard: max. 24 / Expanded: max. 32.
Volume range	Varispan: 0.5 µL – 5 mL Main Array: 1- 200 µL for 96-channel; 0.5 – 25 µl for 384-channel	Varispan: 0.5 µL – 5 mL MDT: Dynamic volume range from 2nL (PinTool) to 235 µL (96Tips P235)
Tip management	Sunken Tip Storage: Recessed deck allows tip boxes to be stored below the work surface.	Surface Storage: Standard deck-level tip placement.
Specialized versions	Fontus NGS: Optimized for the latest genomics workflows. Fontus flow cytometry cocktail prep workstation	JANUS BioTx: Dedicated for automated protein purification/chromatography.
Lab footprint	Efficient: The Standard platform (54") provides high capacity in a streamlined frame.	Scalable: The Expanded platform (68") is for those who need the absolute largest workspace

Which model should you choose?

The JANUS G3 systems support workflows requiring specialized modules or integration with an external device without the addition of an external robotic arm. If your lab's primary focus, for instance, is Protein Purification, the JANUS BioTx workstation provides an "out-of-the-box" solution.

The Fontus systems support workflows requiring throughput and deck density. If you are running high-volume molecular biology or cell-based screening, the Fontus workstation's ability to store more tips and plates in a smaller footprint combined with the Main Array Gripper will significantly reduce your total "time-to-result."

Finding your ideal instrument

While technical specifications provide a roadmap, we understand that every laboratory's workflow has its own unique nuances. Whether you are navigating complex reagent viscosities, strict clean-room requirements, or the need for future-proof scalability, our application specialists are here to help you weigh these variables.

We invite you to contact your local Revvity representative for a personalized workflow consultation. By discussing your specific volume ranges, labware types, and throughput goals, we can assist you configure the exact workstation that will help accelerate your research today and adapt to your discoveries tomorrow

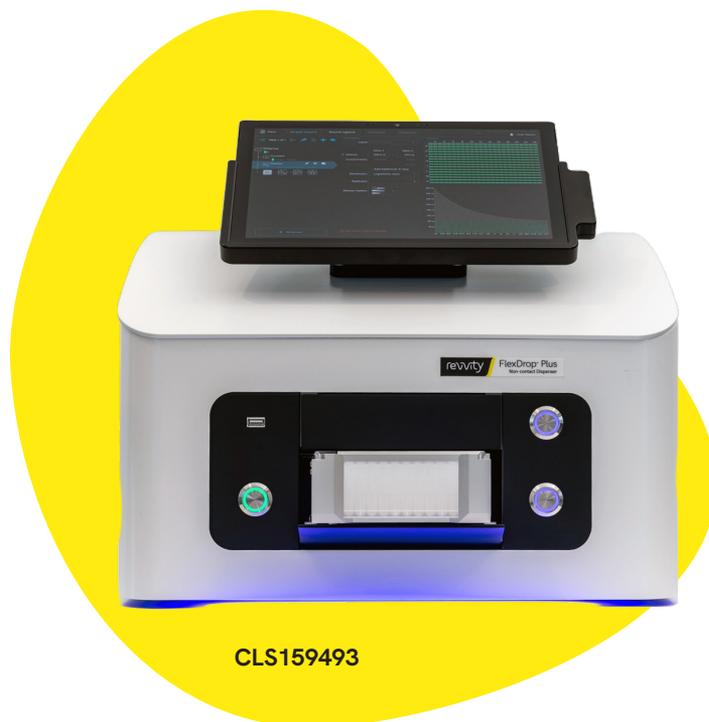
FlexDrop Plus dispenser – contactless nanoliter precision

The FlexDrop™ Plus dispenser enhances your lab's efficiency and budget with its dispensing technology that delivers research-grade precision while helping to reduce reagent costs and workflow bottlenecks.

- Wide volume range (8 nL-80 µL) allows compounds & buffer backfilling on the same instrument
- Low dead volume <1 µL
- Make, optimize and dispense a wide range of liquid classes, incl DMSO
- Wide range of sample types, large and small molecules
- Non-contact positive pressure technology supports the reduction of process cross-contamination
- Fast, reliable, and simple – dispense 96 wells in 10 seconds
- Precision less than 5%CV at 50 nL volume (H₂O) with built-in droplet verification for reliable dispense

The FlexDrop Plus dispenser is especially suited for these applications:

- **Assay development** – Precise gradient dispensing for ELISA, dose-response curves
- **High-throughput screening** – Complex drug concentration matrices and DoE
- **Genomics and proteomics** – Low dead volume for precious samples
- **NGS library prep and qPCR** – Reaction miniaturization and normalization/pooling



Key benefits

- **Cost savings through miniaturization** – Reduce reagent usage by up to 10x, especially valuable for expensive assays
- **Speed and efficiency** – 8-channel simultaneous dispensing with intuitive touchscreen programming
- **Accuracy and precision** – CVs < 5.0% for volumes >100 nL with physical droplet verification
- **Versatility** – Handles aqueous, DMSO, glycerol, beads, and high-viscosity solutions across 96/384/1536 well plates

Resources:

- [FlexDrop Plus Dispenser specifications sheet](#)
- [Revolutionizing ADC development: The critical role of advanced dispensing systems](#)
- [FlexDrop Plus overview video](#)



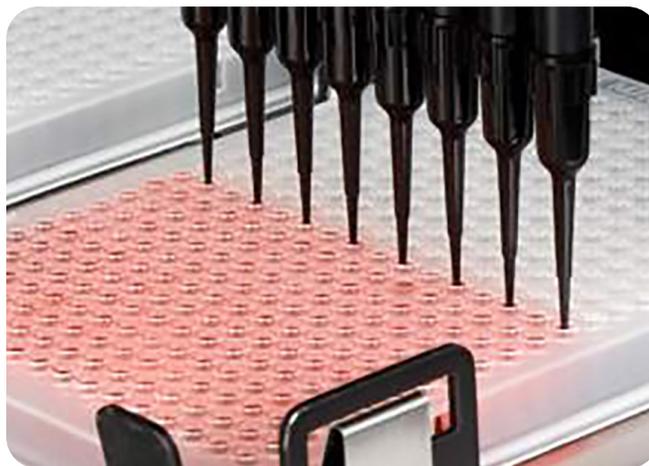
Learn more

More information
and resources

SECTION 2

The science of the dispense

Revvity offers three pipetting technologies to provide capabilities to address the challenges posed by the physical properties of the reagents used in protocols.



Revvity pipetting technology decision matrix

Technology	How it works	Recommended for	Key Advantage(s)	Primary instrument(s)
Air displacement	Uses an air cushion between the internal piston and the liquid	Aqueous liquids, DNA/RNA prep, standard buffers.	Disposable tips prevent carry-over	<ul style="list-style-type: none"> AssayMate JANUS with MDT head Fontus with 96/384 head Zephyr
Liquid displacement	Piston acts on a system liquid, eliminating the air cushion.	Viscous (oils/detergents), Volatile, or foaming liquids.	Improved reproducibility for "difficult" liquids that cling to standard tips.	<ul style="list-style-type: none"> Fontus and JANUS G3 with Varispan head
Non-contact	Reagents are "jetted" via pressure without tip contact.	Reagent conservation, HTS spiking, sensitive biologics (ADCs).	Nanoliter precision reduces reagent costs (<1 μ L dead volume); non-contact prevents cross-contamination; eliminates tip costs.	<ul style="list-style-type: none"> FlexDrop Plus
Mixed / context	Combines multiple technologies on a single deck.	Complex, multi-stage workflows (e.g., Protein Purification).	High flexibility enabling the use of the optimal technology for each reagent in one run.	<ul style="list-style-type: none"> explorer G3 JANUS BioTx

Please [contact us](#) to discuss your own requirements.

SECTION 3

Applications, the solution map

Revvity’s liquid handling portfolio is not a one-size-fits-all solution. We map our technology to your specific scientific milestones, ensuring that your instrument choice supports your current research while providing a pathway for future expansion.

For NGS solutions, please consult our [NGS dedicated brochure](#).



Guide by application

Application Type	Primary Task	Throughput Need	Primary Instrument (s)
Routine Lab Tasks	Serial dilutions, ELISA, Plate Coating/Washing	Low to Mid	AssayMate workstation
Molecular Biology	DNA/RNA Extraction, PCR setup, Sequencing Prep	Mid to High	Zephyr G3, Fontus, JANUS
Cell Biology	Cell-based assays, Organoid research, Seeding	Low to Mid	AssayMate, Fontus, JANUS
Cytometry / Staining	Antibody cocktailing, Cell surface staining	Mid to High	Fontus cocktail prep workstation FlexDrop Plus
Protein Sciences	Crystallization Screening, Purification, Immunoprecipitation (IP)	Low to Mid	Fontus, JANUS
Biologic Screening	ADCs, Nanoliter “Spiking”, Dose Response	High	FlexDrop Plus
Drug Discovery	Primary Reformatting, Lead Optimization, ADME	High (Versatile)	Fontus, JANUS
Solid Phase Extraction	Toxicology, Food/Soil analysis, Mass Spec Prep	Mid to High	Zephyr G3 SPE
Industrial Screening	Integrated HTS, 24/7 Automated Workstations	Ultra-High	explorer G3

Please [contact us](#) for more information.

Guide by sample type

Sample Type	Primary Instruments	Key Considerations
Small Molecules	FlexDrop Plus, Fontus, JANUS	DMSO compatibility, acoustic dispensing, high throughput
ADCs	FlexDrop Plus	Nano-to-microliter dispensing, DMSO compatibility, contamination prevention, non-contact dispensing to preserve conjugate integrity
Biologics	FlexDrop Plus, Fontus, Zephyr, JANUS	Temperature control, gentle handling, viscous liquid capability
Cells & Tissues	AssayMate, Fontus, JANUS, Zephyr	Sterile technique, gentle dispensing, CO ₂ incubator integration
Nucleic Acids	AssayMate, Fontus, JANUS, Zephyr	Contamination prevention, magnetic bead handling, precision
Complex Matrices	Zephyr G3 SPE, JANUS G3 SPE	Solvent resistance, regulatory compliance, audit trail

Please [contact us](#) for more information.

Choosing your ideal instrument

Identifying your application requirements is only the first step in the journey toward optimized laboratory automation. While the table above serves as a foundational guide for matching scientific tasks with Revvity's specialized hardware, the true power of our portfolio lies in its customizability. Whether you are automating Dose Response studies on a FlexDrop Plus dispenser or performing complex

Solid Phase Extraction on a Zephyr G3 workstation, the ideal configuration depends on your specific labware, your required dynamic range, and your daily throughput goals.

Because no two workflows are identical, even within the same application, we invite you to leverage our deep technical expertise. Contact us today for a personalized workflow assessment.

Resources:

Visit our [Resource Center](#) to browse the resources for the liquid handling applications.



Learn more

More information
and resources

Application-dedicated workstations

Additionally to our general liquid handling portfolio, application specific workstations have been configured to help address specific workflow requirements of contemporary drug discovery and diagnostics research, including biotherapeutics, genomics and drugs of abuse detection.

For NGS dedicated workstations, please consult our [NGS dedicated brochure](#).

Workstations	Ref.	Dedicated application	Main advantage
Zephyr™ G3 SPE	CLS145303 , CLS145304	Solid Phase Extraction	Clog Detection: Integrated ultrasonic sensors identify clogged wells in real-time to preserve precious samples. It helps prevent the risk of clogged sample wells overflowing into next sample wells, thus enabling full walkaway.
Fontus™ cytometry cocktail prep	NTUSCOCKTAILPREP	Flow Cytometry Prep	Standardization: Automates antibody cocktailing and staining with digital audit trails for total traceability.
JANUS™ G3 BioTx	YJS80P1 , YJM8MP1	Protein Characterization	Multi-mode Purification: Supports column, tip, and batch chromatography on a single platform to accelerate process development.

Resources:

- [Fontus flow cytometry cocktail preparation flyer](#)
- [Fontus webinar: precision and efficiency in flow cytometry workflows](#)



Learn more

More information and resources

REVVITY AND BIOLEGEND ON-DEMAND WEBINAR:

Unlock precision and efficiency in your flow cytometry workflows.

Watch now

SECTION 4

Better together – platform synergy

Bulk preparation and nanoliter precision

Pair the high-volume capacity of a Fontus, JANUS, or Zephyr liquid handler with the nanoliter precision of a FlexDrop™ Plus dispenser. This synergy allows for bulk master-mix preparation followed by ultra-low volume “spiking,” potentially up to 90% in reagent costs.

In a typical high-throughput workflow, reagents fall into two categories: high-volume buffers/diluents and high-value triggers/samples. By pairing a Revvity liquid handler with the FlexDrop™ Plus dispenser, you can use the right tool for each task to help optimize your budget and your data quality.

1. Potential drastic reagent savings through miniaturization

The FlexDrop Plus enables “tipless” dispensing with a dead volume of less than 1 µL. While a standard liquid handler prepares your master mixes or reformats primary samples, the FlexDrop Plus can dispense expensive enzymes, antibodies, or compounds in volumes as low as 8 nL.

- **The Benefit:** Transition from 20 µL reactions to 2 µL (or less). Many labs report reducing reagent consumption by up to 90%, allowing you to run 10x more samples with the same kit budget.

2. Helps prevent cross-contamination with “Tipless” technology

The FlexDrop Plus uses positive pressure to “jet” droplets directly into the destination plate without any physical contact.

- **The Benefit:** Eliminate the recurring cost of pipette tips for your most frequent dispenses. This contactless approach is ideal for sensitive biologics (ADCs) and PCR master mixes where purity is non-negotiable.

3. Overcoming physical fluid limitations

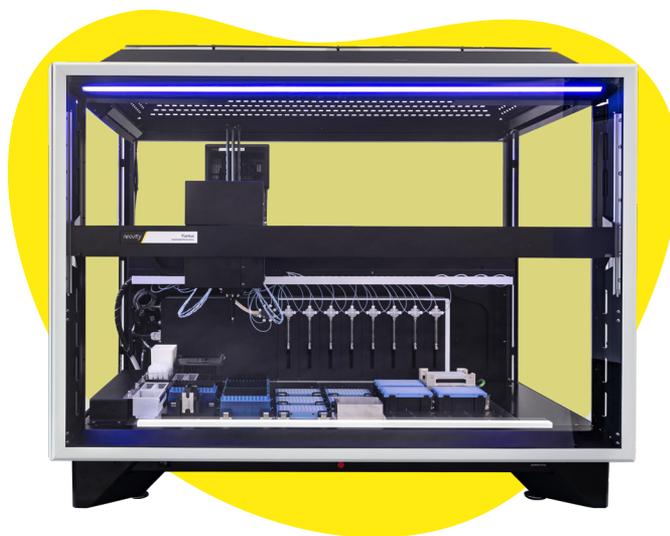
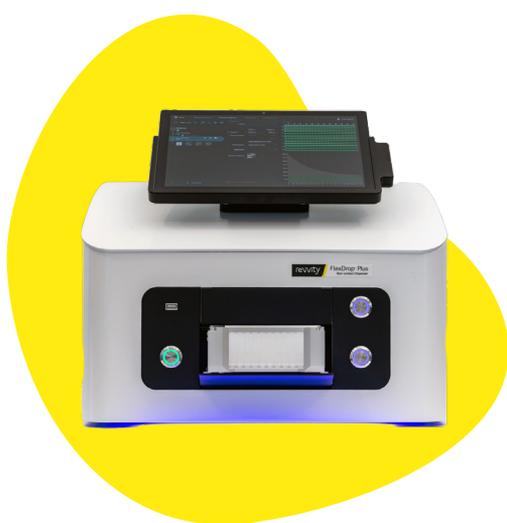
Revvity liquid handlers (JANUS/Fontus) excel at complex movements and viscous liquids via Liquid Displacement. The FlexDrop Plus complements this by handling a vast range of liquid classes (Aqueous, DMSO, and up to 50% Glycerol) with integrated droplet verification.

- **The Benefit:** The FlexDrop Plus uses built-in sensors to verify that droplets are dispensed as intended. If a well is missed, the system can automatically re-dispense, supporting data integrity for your most critical assay triggers.

4. Application Spotlight

Application	Revvity Liquid Handler Role	FlexDrop™ Plus Role
NGS Library Prep	Bead clean-up and pooling	Low-volume enzyme/adaptor addition
qPCR Setup	DNA/RNA template normalization	Master mix and primer dispensing
Drug Discovery	Primary sample reformatting	Nanoliter dose-response “spiking”
Cell Assays	Media exchange and cell seeding	Compound or cytokine stimulation

Don't let high reagent costs or tip consumption limit the scale of your research. By integrating the FlexDrop™ Plus into your automation strategy, you can gain the ability to scale down volumes enabling you to scale up your discovery potential.



Resources:

Application available soon

Plug-and-Play Miniaturized Electroporation: A Seamless Integration of the DropGenie Platform into the Revvity Fontus and FlexDrop Liquid Handling Ecosystems for High-Efficiency Cellular Delivery

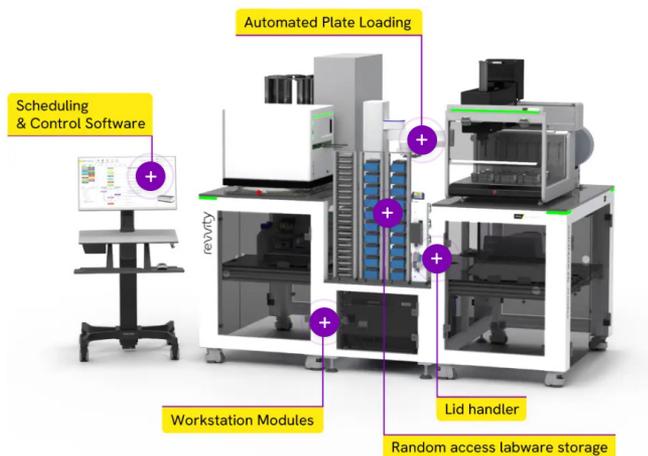
SECTION 5

explorer G3 workstation: Integrated automated solutions

Full workflow automation

From cellular screening and imaging applications to high-throughput screening and genomics-based applications, custom explorer™ G3 integrated workstations provide innovative application-focused laboratory automation solutions which simplify microplate handling, liquid handling, and detection.

- **The power of integration:** Scalable solutions ranging from simple plate feeding to complex, multi-level explorer™ G3 workstations.
- **Intelligent scheduling:** plate::works™ and job::manager™ provide real-time decision-making and a calendar-style view for multi-user environments.
- **Scientific depth:** Application-focused integration for HTS, automated ELISA, and high-content cellular screening and many more.



Custom workstations - up to 8 instruments

[Check this interactive infographic](#)

With complete high-throughput multimode plate detection, high content analysis, and liquid handling portfolios, we are positioned to understand your scientific and application needs. You can count on us to provide solutions that help address the full breadth of your scientific requirements, all from a single source who can deliver the application support you need. Additionally, a variety of instrumentation from other suppliers can be integrated into an explorer G3 integrated workstation.

Please [contact us](#) to discuss your own requirements.

Spectrum of solutions:



Plate feeding - up to 2 instruments
The SmartBench workstation



Workstations - up to 4 instruments



Custom workstations - up to 8 instruments



Custom platforms - up to 20 instruments

Resources:

- [Embracing modularity: flexible automation's role in drug discovery research](#)
- [Partnering with purpose: Australian Organoid Facility From hits to leads: automating lead optimization - SAR assay development scalable solutions](#)



[Learn more](#)

**More information
and resources**

SECTION 6

Why Revvity? Your partner in discovery.

Full workflow automation

From your first benchtop workstation to a fully automated integrated robotic solution, Revvity provides an array of paths for growth tailored to your specific needs. Our modular hardware and application-specific expertise aim to support your laboratory at the forefront of discovery, regardless of your throughput or complexity.

Revvity's commitment is to help transform your scientific challenges into opportunities—from cell and gene therapy to small molecule discovery. We provide a streamlined ecosystem that guides researchers through each critical stage, addressing current obstacles while anticipating the emerging challenges of tomorrow.

1. Platforms and automation: Enhanced success rates

Unlock access to advanced technology designed for reliability. Our instrumentation doesn't just move liquid; it helps move your science forward:

- **Scalability:** Meet a wide array of laboratory needs, from compact benchtop units to high-throughput integrated workstations.
- **Efficiency:** Minimize time-to-market by automating manual steps, integrating data, and streamlining compliance.
- **Precision:** Drastically reduce handling errors through robust readouts and advanced integrated analytics.
- **Innovation:** Tackle unsolved research challenges with validated QC methods for novel therapeutic modalities.



Learn more

Automated
instrumentation

2. Reagents: The fuel for your results

An automated system is only as good as the chemistry it runs. Revvity offers a wide-ranging portfolio of reagents optimized for integration with our platforms:

- **Comprehensive selection:** Immunoassays (AlphaLISA™, HTRF™), gene editing tools (Dharmacon™), and high-quality antibodies (BioLegend).
- **Workflow transformation:** Specialized molecular biology solutions and cell-painting kits that help turn raw samples into actionable data.



Learn more

Discover our
reagent portfolio

3. Revvity Signals: Turning data into insight

Bridge the gap between the wet lab and the digital world. Revvity Signals software solutions serve scientists across Pharma, Biotech, and Clinical Sciences. They integrate generative AI and advanced natural language processing to streamline user workflows.

- **Signals Notebook One™** is a SaaS Electronic Laboratory Notebook that natively integrates ChemDraw™ and Spotfire®, an end-to-end R&D workflow solution providing data capture and collaboration, data processing, and data-driven analytics. Signals One sets the stage for AI success by putting customers' data in context with AI-ready data management.
- **Signals VitroVivo:** Streamline the analysis of complex assays with automated calculation templates and robust data integration.



Learn more

Learn more.

4. Revolutionary support: Maximizing your uptime

We believe your success is entwined with ours. Revvity provides more than just a repair service; we provide a proactive collaboration to keep your lab running at peak efficiency.

- **Expert response:** Our Field Service Engineers and technical support teams are renowned for rapid response times and high first-time fix rates.
- **Specialized services:**
 - **IQ/OQ:** Comprehensive instrument qualification for regulated environments.
 - **Adept support:** On-demand application and software training tailored to your team.
 - **Proactive care:** Scheduled preventative maintenance to eliminate downtime before it happens.



[Learn more](#)

Review our maintenance services.

Challenge accepted.

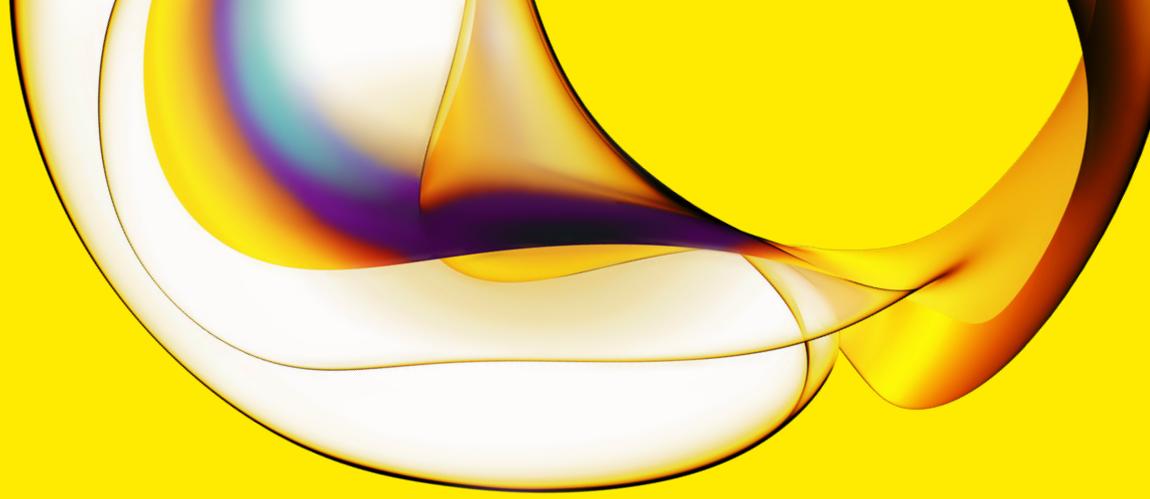
At Revvity, we don't just provide tools; we provide the bridge between a question and a breakthrough. Whether you are looking to automate a single assay or build the lab of the future, we are ready to help you navigate the path to impactful discovery.

Welcome to the future of science.
Welcome to Revvity.



[Learn more](#)

Contact an Expert



www.revivity.com

revvity

Copyright ©2026, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

2538600

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