



Achieve peak performance in your genomic workflow

At Revvity, our optimized, automated solutions are designed to improve the efficiency of your genomic workflows. By providing labs with complete, single-source solutions from sample to analysis, we aim to address the complex challenges associated with genomic analysis.

Custom solutions that work for you

Sample disaggregation
Omni homogenizers



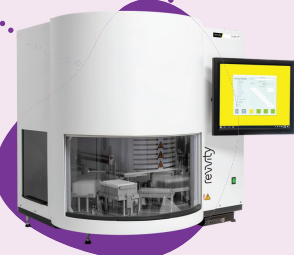
Quantitation of nucleic acids
Plate reader



Library prep
Automated liquid handling



Library prep kits



Nucleic acid isolation
chemagic™ technology



Visual QC & sizing of nucleic acids
Microfluidic devices



Analysis
Microbiome analysis

Sample homogenization

Bead mill homogenizers

The Omni Bead Ruptor bead mill homogenizers leverage the power of mechanical bead-beating to maximize tissue homogenization and cell lysis, with high-impact processing speeds up to 8 m/s.



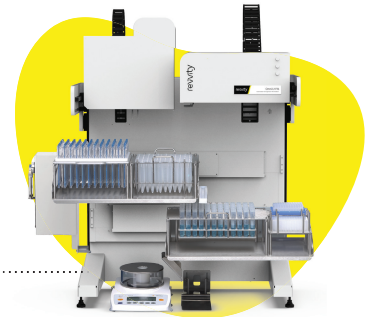
Rotor stator homogenizers

Handheld or stand-mounted high-shear laboratory homogenizers designed to process tough samples quickly and efficiently without the need for harsh chemical lysis. Together with faster homogenization times that minimize heat degradation, when coupled with single-use OmniTip™ disposable probes that significantly reduce risk of contamination, sample integrity is maintained.



Automated homogenizer

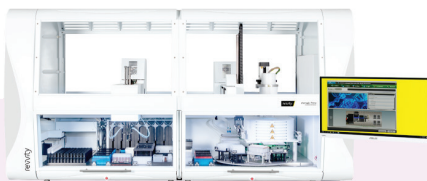
Omni automated homogenizer workstations enable labs to future-proof their throughput with not only robotic homogenization, but the entire front-end sample prep for tissue lysis, including sample weighing, buffer addition/liquid handling, and tube-to-plate reformatting. Be confident automating your workflows from the first step.



Nucleic acid isolation

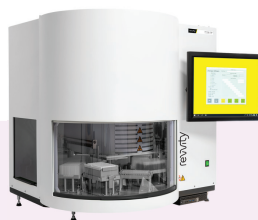
The chemagic™ instruments are flexible and automated systems for high-quality DNA/RNA isolation based on proven M-PVA Magnetic Bead Technology.

- The magnetized rods transfer magnetic beads through various process solutions without transferring liquid, effectively minimizing the risk of sample cross-contamination.
- Rotation of the demagnetized rod heads efficiently resuspends the particles during washing steps, enabling high yields, purity, and low levels of fragmentation of nucleic acids.



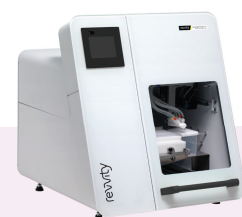
chemagic Prime

- Sample volumes from 50 µL - 10 ml
- 1-192 samples per run
- Tracked reagent/tip consumption and UV disinfection
- On-deck heating/cooling/shaking options



chemagic 360

- Sample volumes from 50 µl - 18 ml
- 12-96 diverse sample materials per run
- 1D/2D barcode scanning



chemagic Prepito™

- Sample volume up to 1 ml
- Up to 12 diverse sample materials per run
- Barcode reading for sample tracking

Microfluidic analysis

If you need to analyze proteins, nucleic acids, or both our LabChip™ automated microfluidic capillary electrophoresis (micro-CE) technology enables the simplification of traditional gel separations, resulting in robust analysis in a fraction of the time.

- Separate, identify and analyze genomic and protein samples in seconds
- Visualize your data as an electropherogram, virtual gel, or tabular report



Library prep kits

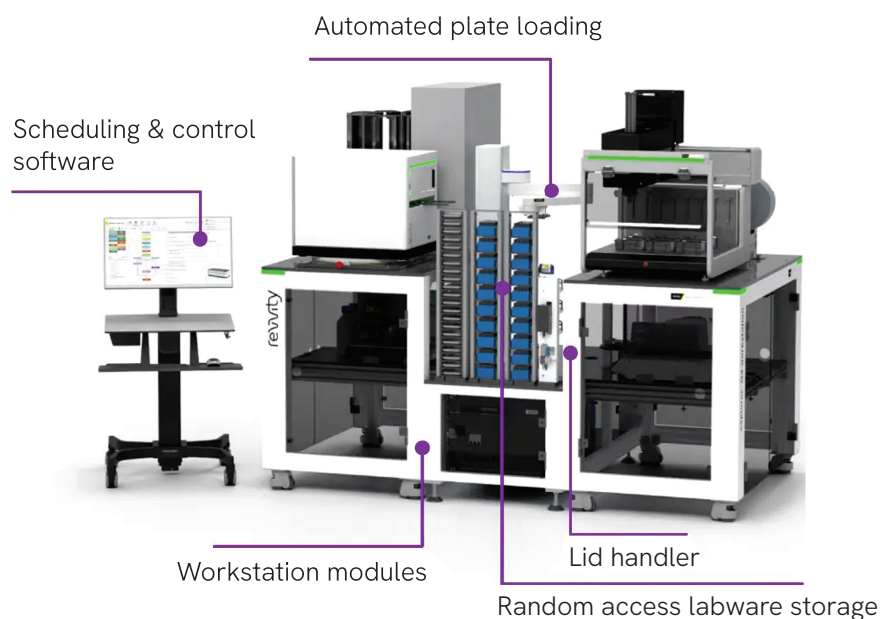
Revvity offers a complete portfolio of NGS library preparation kits and barcodes designed to achieve robust results every time, from low to high-throughput applications. The NEXTFLEX™ kits cover most applications such as whole genome sequencing, gene expression analysis, small RNA, metagenomics and agrigenomics.



Integrated laboratory automation

Custom explorer™ G3 integrated workstations provide innovative application-focused laboratory automation solutions which simplify microplate handling, liquid handling, and detection. Additionally, a variety of instrumentation from other suppliers can be integrated and custom platforms up to 20 instruments

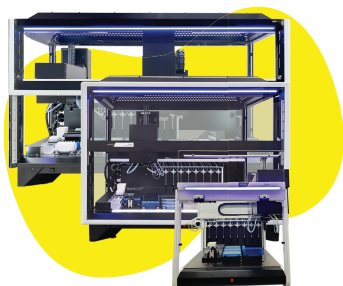
- Automated plate loading
- Automating cellular, drug discovery, and genomics workflows
- Workstation modules
- Scheduling and control software
- Robotics



Automated liquid handling workstations

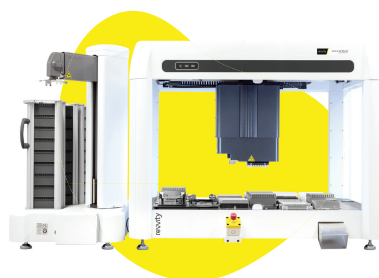
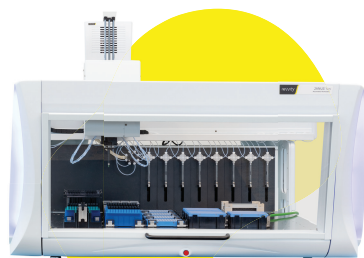
Fontus

With easy-to-use software, optimized deck access, verified protocols, and reformatting capabilities to simplify your workflows and improve turnaround times, the new Fontus™ liquid handler is a powerhouse, helping you to achieve even more.



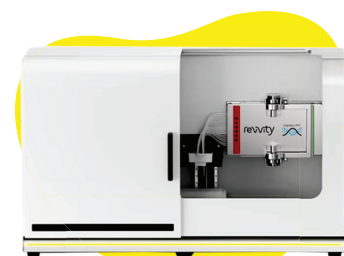
JANUS

The JANUS™ G3 workstation can be configured based on the size of the instrument, pipetting arm technology, labware movement options, and a number of other accessories to fully automate your application. To accommodate your choice of deck size and pipetting arm, Revvity offers four different versions of the JANUS automated workstation.



Sciclone

With proven performance, open-deck design, and integration-friendly architecture, Revvity's Sciclone™ G3 liquid handling workstations offer 80+ vendor-qualified automated NGS methods to enhance your workflow efficiency.

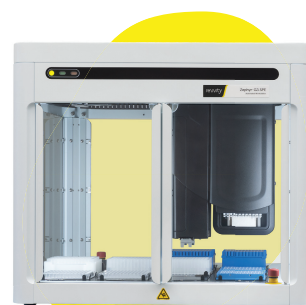


BioQule NGS solution

The BioQule™ NGS solution automates DNA library preparation, supporting up to 8 samples per run with minimal manual intervention. Compatible with leading library prep chemistries, it integrates DNA extraction, library prep, and quantification into a streamlined workflow. With an intuitive interface and up to 80% reduced hands-on time, it delivers robust, high-quality libraries for Illumina sequencing, making it ideal for low-throughput labs.

Zephyr

The Zephyr™ G3 workstation's small footprint makes it ideal for workbench operation, while the convenient deck design provides ready access from all four sides to consumables and accessories. The Zephyr G3 workstations offer 40+ automated vendor-qualified NGS methods.



FlexDrop Plus non-contact dispenser

The FlexDrop™ Plus system dispenses a wide range of reagent and sample types fast with high accuracy and low dead volume. Small bench footprint too!



revvity