



Accelerate your time to market with OEM and customization services.

Why partner with Revvity?

We don't believe in one-size-fits-all. Every OEM customer receives dedicated, one-on-one support tailored to their unique needs – because your challenges deserve personalized attention, not a generic response.

Deep scientific expertise

Decades of genomic workflow knowledge

Personal, dedicated OEM support

A single point of contact who knows your project inside and out

End-to-end partnership

From R&D and feasibility studies to serial manufacturing and delivery

Trusted globally

Supporting genomic applications worldwide

Custom formulation and fill volumes

We customize kit formats to match your workflow and product requirements.

Customized private labeling and branding

From volumes to packaging formats, we adapt to your production and distribution needs.

Custom kit configurations

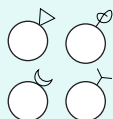
Your brand, your identity – we support white-label and co-branded labelling solutions.



Your trusted partner from sample to sequence.

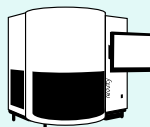
Sample preparation

Magnetic beads



Our polyvinyl alcohol based magnetic beads can be functionalized to bind your specific target molecules, providing the foundation for reliable, consistent results across diverse applications.

Nucleic acid purification platforms



Flexibility is built in. Our chemagic™ platforms seamlessly integrate into your workflow and scale with your business, whatever your throughput demands today or tomorrow - the platforms cover both a wide range of sample volume (50 μ L to 18 mL) and different levels of throughput (12/24/96) with high-quality DNA/RNA yields.

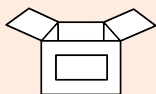
Nucleic acid purification chemistry



Our extensive portfolio offers reliable DNA and RNA isolation solutions for diverse sample types. In today's rapidly evolving genomic research landscape, we continuously optimize our chemistry to adapt to emerging applications.

NGS library preparation

Library prep kits



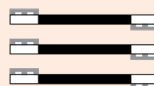
DNA and RNA kits enable whole-genome sequencing, mRNA, total RNA, strand-specific sequencing, small RNA, and microRNA analysis, empowering researchers to study genetic variations, gene expression, and regulatory mechanisms.

Barcodes



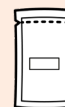
Revvity offers high-performance NGS solutions, including barcoded adapters with unique dual indexes (UDI) compatible with Illumina®, Watchmaker™, Twist®, and other TruSeq®-style kits, plus adapters combining UDIs with unique molecular identifiers (UMI).

Universal blockers



Our universal blockers enhance targeted sequencing workflows, delivering market-leading performance at a fraction of the cost of major competitors.

Accessories



Our accessory products, include NGS Cleanup Beads, poly(A) Enrichment Kits, and ribodepletion tools are compatible with most major library prep protocols, offering cost-effective and flexible solutions to enhance workflows while maintaining performance.

Ready to bring your vision to life? Revvity's OEM expertise, proven technology, and end-to-end flexibility make us the partner you've been looking for.

Let's build something together.



Learn more

Scan or click
for more information

revvity