



RNA-seq solutions portfolio.

Comprehensive tools for every RNA-seq need

RNA sequencing (RNA-seq) is the gold standard for capturing the dynamic expression landscape of cells. Revvity's comprehensive portfolio—from library preparation kits to advanced enrichment, depletion, and multiplexing solutions—streamlines your workflow and enhances the detection of low-abundance transcripts.

Our comprehensive portfolio includes:

- **Complete workflow integration**
Seamlessly designed for use with Illumina® and Element® sequencing platforms, our modular products ensure directional library preparation and accurate transcript mapping.
- **Enhanced mRNA purification**
NEXTFLEX™ Poly(A) Beads 2.0 use magnetic bead-based technology for high mRNA yield, minimal rRNA carry-over, and compatibility with both manual and automated systems.
- **Efficient ribodepletion solutions**
Our ribodepletion kits, available in CRISPR-based or hybridization-probe formats, remove >80% of rRNA to improve the detection of rare, biologically relevant transcripts.
- **Robust multiplexing with advanced barcoding**
NEXTFLEX RNA-seq 2.0 UDI Barcodes and NEXTFLEX UDI UMI Barcodes offer error correction, prevent index mis-assignment, and incorporate Unique Molecular Identifiers (UMIs) for precise quantification.

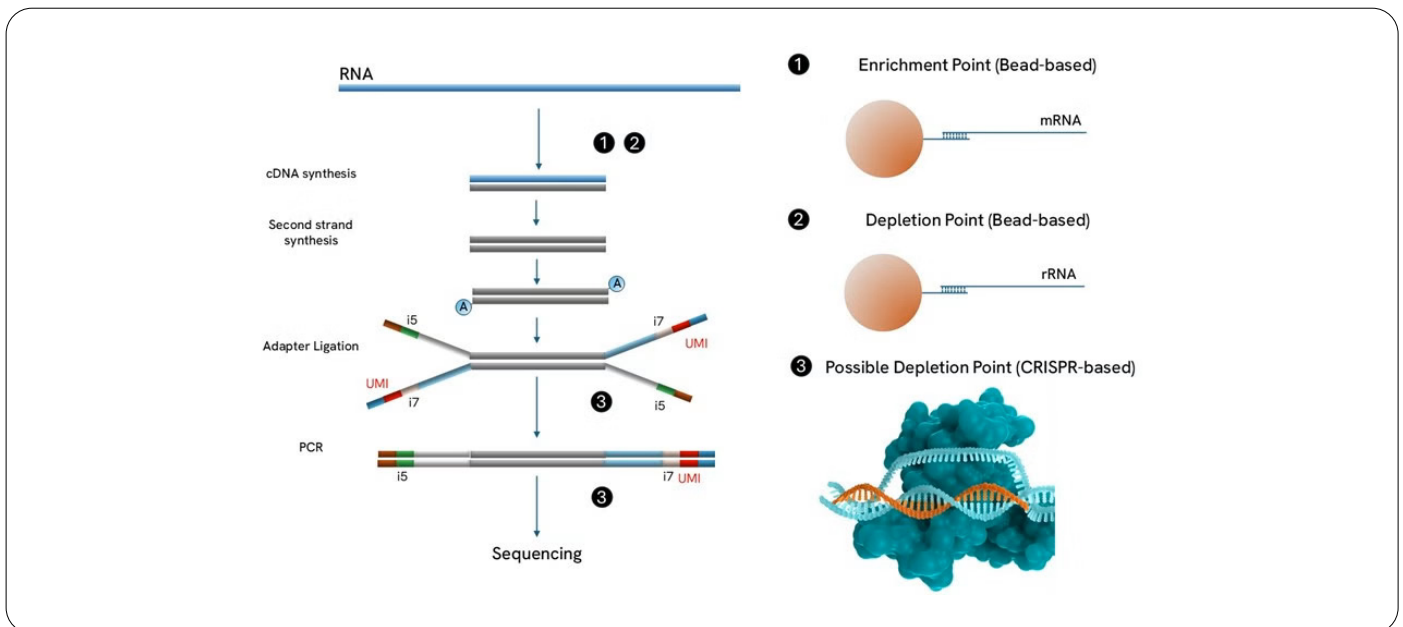


Figure 1: This figure illustrates the classic RNA-seq workflow. The numbers in the diagram show where various enrichment & depletion based accessory products can be used in RNA-seq workflows.

RNA-seq library preparation and mRNA enrichment

NEXTFLEX RNA-seq library prep kits

- Convert RNA into high-quality, directional cDNA using optimized workflows.
- Preserve strand information for accurate transcript mapping.

Product Spotlight: NEXTFLEX Poly(A) Beads 2.0

- **Key features:**
 - Delivers improved mRNA yield with low rRNA contamination
 - Magnetic bead-based isolation for 10 ng – 5 µg of total RNA.
 - Suitable for both manual and automated isolation.

Multiplexing with advanced barcoding

NEXTFLEX RNA-seq 2.0 UDI Barcodes

- **Highlights:**
 - 384 unique dual-index adapters engineered for robust multiplexing.
 - Minimizes index hopping and mis-assignment on patterned flow cells.

NEXTFLEX UDI-UMI Barcodes

- **Highlights:**
 - Integrate Unique Molecular Identifiers (UMIs) to distinguish true PCR duplicates from genuine molecules.
 - Improve quantitation, variant calling, and overall library complexity assessment.

Advanced ribodepletion solutions

Our approach:

- Choose from CRISPR-Cas9-based or hybridization-probe-based kits to remove rRNA and unwanted transcripts.
- **Benefits:**
 - Increased detection of rare transcripts
 - Reduced background noise
 - Enhanced statistical power for gene expression studies

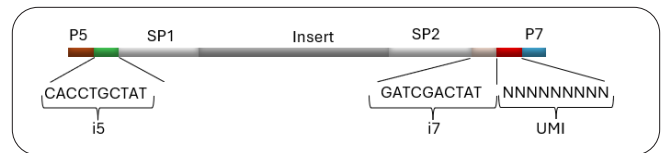


Figure 2: A diagram showing UMI layout on a library molecule.

Transform your RNA-seq experiments today.



LEARN MORE

revvity