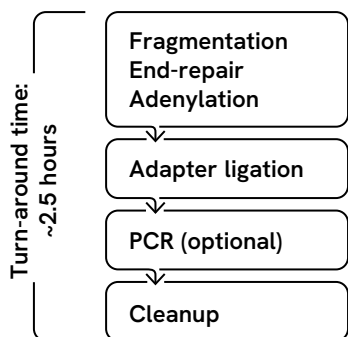




High-performance library prep for standard and miniaturized workflows.

Streamlined library prep

The new NEXTFLEX™ Rapid XP v3 DNA-seq kit enables library preparation from a broad range of sample types, including low-input and FFPE DNA samples. This kit is optimized to convert high- and low-quality DNA into Illumina® and Element® compatible libraries with improved yield, low sequencing artifacts and uniform coverage. It can deliver PCR-free libraries in under 90 minutes and scales easily to high-throughput workflows and automation. A miniaturization protocol is available for price-sensitive applications.

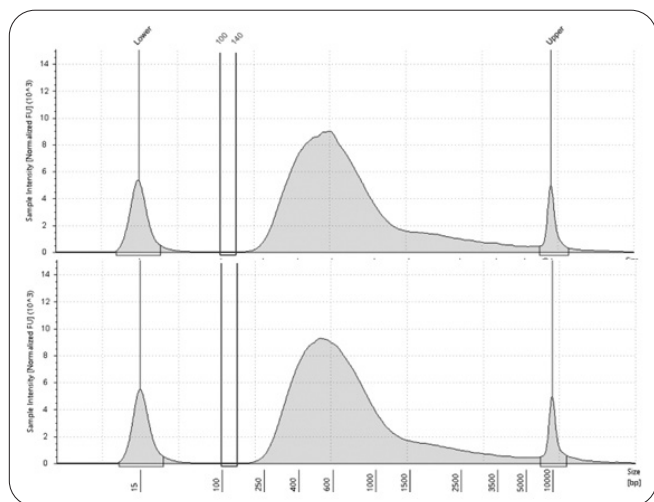


NEXTFLEX Rapid XP v3

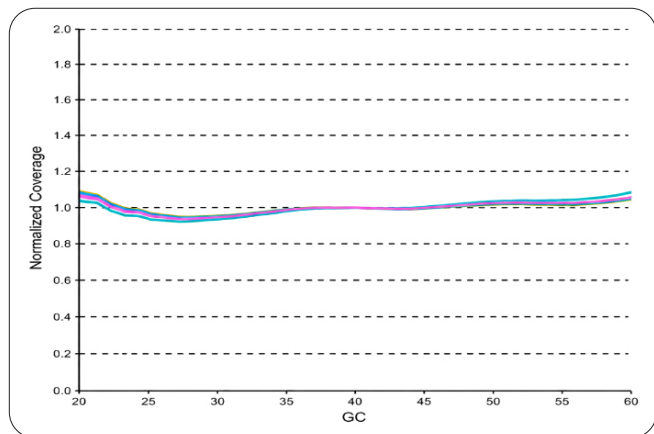
- Working range from <1 ng – 500 ng.
- Enzymatic fragmentation with end-repair and A-tailing in one reaction
- Highly efficient ligation and uniform coverage
- PCR-free workflow with DNA inputs ≥ 75 ng.
- Up to 1,536 UDI barcodes available for scaling and high-throughput multiplexing.
- Automated on the Sciclone™ G3 NGSx, Sciclone G3 NGSx iQ™ and Fontus™ NGS workstations.
- Miniaturization protocol available for the FlexDrop™ Plus Dispenser (Revvity) and I.DOT Liquid Handler (Dispendix®)
- Compatible with NEXTFLEX™ Normalization Beads to shorten time needed for quantification and pooling of final libraries.

Robust performance even at miniaturized volumes

The NEXTFLEX Rapid XP v3 DNA-seq kit offers a streamlined workflow with highly efficient ligation, generating low adapter-dimer levels and consistent library yields even at miniaturized reaction volumes.



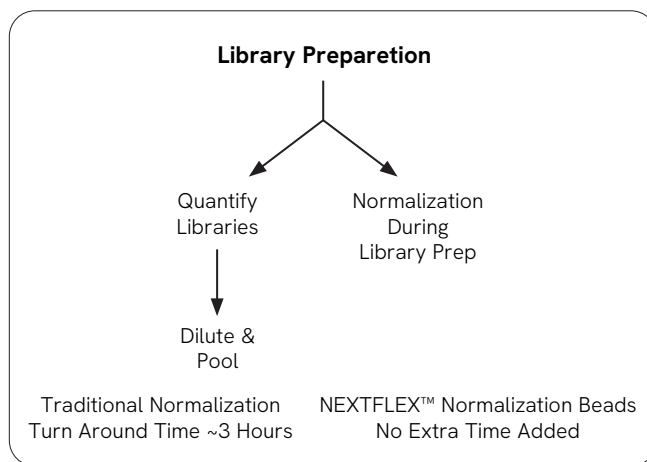
Consistent yield and low dimer formation with NEXTFLEX Rapid XP v3 DNA-seq kit. Different library replicates were prepared using 25 ng of human gDNA as input, at one-quarter volume reaction. Adapter dimer should appear as a peak at ~140 bp.



Analysis of GC bias from replicate libraries obtained with NEXTFLEX Rapid XP v3 DNA-Seq kit. Results indicate a consistent coverage across 20 to 60% GC content.

Bead-based normalization option

Libraries prepared with the NEXTFLEX Rapid XP v3 are compatible with NEXTFLEX™ Normalization Beads. Beads are sold separately.



Compatibility with target enrichment workflows

Libraries prepared with the NEXTFLEX Rapid XP v3 DNA-seq kit are compatible with downstream hybridization-based target enrichment workflows. This provides flexibility to use probes from multiple vendors, supporting applications such as custom targeted sequencing and exome sequencing.

Flexible multiplexing options for high-throughput applications

We offer a range of barcodes compatible with the NEXTFLEX Rapid XP v3 DNA-seq kit to improve multiplexing capabilities for both low and high-throughput multiplexing needs. A set of up to 1,536 NEXTFLEX UDI adapters are now available. A 96-barcode NEXTFLEX UDI-UMI set is also available for detection of low frequency variants. Barcodes are sold separately. Visit our website for compatible barcode options:



[Learn more](#)

Scan or click to
visit our website

revvity