



Library prep + normalization = faster, easier, library prep at a lower price point

Fully automated normalization and fragmentation

The new NEXTFLEX® Rapid XP V2 DNA-seq kit includes proprietary NEXTFLEX® normalization beads*, which provide a consistent mass and sequencing cluster density for all samples in a library pool. This shortens the time needed for quantification and pooling preparation for sequencing in a high throughput lab by up to 3 hours per 96 samples.

**The kit developed for BioQule™ NGS System does not include normalization beads but the instrument quantifies each library before completion of the workflow*

Fragmentation
End-Repair
Adenylation

Adapter Ligation

Cleanup

Post-Ligation PCR

Normalization

NEXTFLEX® Rapid XP V2 DNA-seq kit for Illumina® platforms

- Fully automated normalization & fragmentation
- Working range from 100 pg - 1 µg
- Enzymatic fragmentation with end-repair and A-tailing in one reaction
- Highly efficient ligation with low adapter dimer formation
- No need to dilute adapters within range of 10 ng - 1 µg
- PCR-free workflow with high DNA inputs
- Efficient: Automated on the Sciclone® G3 NGSx, Sciclone® G3 NGSx iQ™ workstations, Zephyr® NGS workstations for mid-high throughput automation, and now on the BioQule™ NGS System for low throughput labs
- Up to 1,536 UDI Barcodes available for multiplexing

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

Highly efficient ligation with low adapter formation

The NEXTFLEX® Rapid XP V2 DNA-seq kit offers a streamlined workflow with highly efficient ligation, generating low amount of adapter dimers even at low input concentrations.

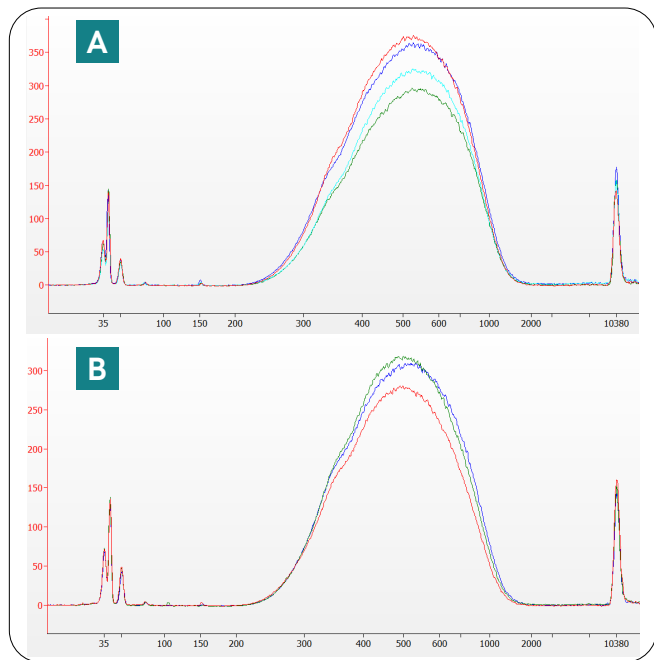


Figure 1. Low dimer formation with NEXTFLEX® Rapid XP V2 DNA-Seq kit. Different library replicates were prepared using either 500 pg (A) or 1 ng (B) of high-quality genomic DNA as input. Adapter dimer should appear as a peak at ~150 bp.

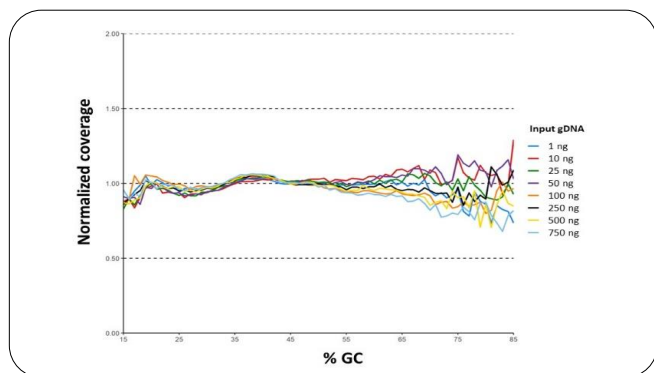
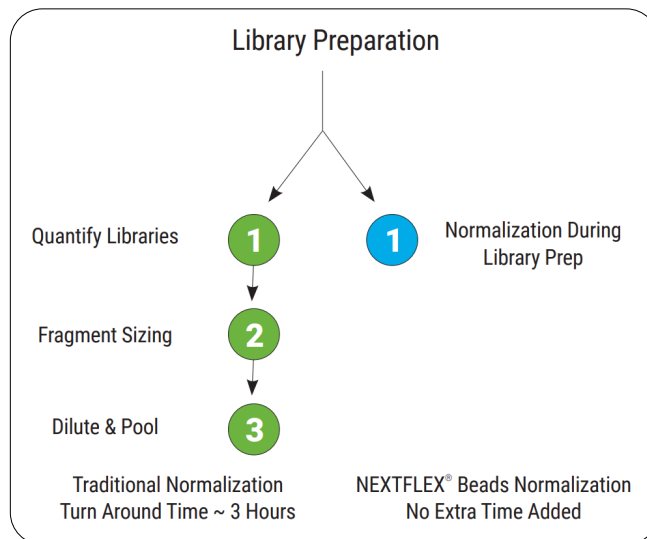


Figure 2. Analysis of GC bias from libraries obtained with NEXTFLEX® Rapid XP V2 DNA-Seq kit. Results indicate a consistent behaviour regardless of input amount from 15 to approximately 60% GC content. Above that some bias is observed, specially at high inputs.

Simplify library construction with NEXTFLEX® normalization beads



Flexible multiplexing options with 1,536 barcodes available for high-throughput multiplexing

We offer a broad range of color-balanced NEXTFLEX® adapters that improve multiplexing capabilities for both low-level and high-level multiplexing needs. A set of 1,536 NEXTFLEX® UDI adapters which are compatible with the NEXTFLEX® Rapid XP DNA-seq kit v2 are now available for ultra high-throughput applications.

Barcodes are not supplied with the NEXTFLEX® Rapid XP DNA-seq kit v2 and need to be purchased separately.

For more information about barcode options compatible with the NEXTFLEX® Rapid XP DNA-seq kit v2 and bundled pricing, please contact www.revivity.com.

Contact an NGS expert and get your DNA project started today!

revvity