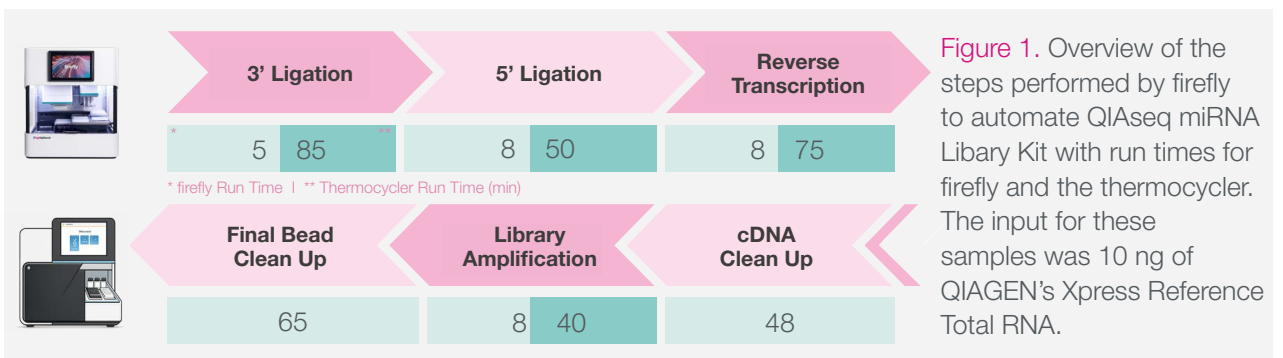


Automating QIAseq[®] miRNA Library Kit on firefly[®]

Introduction

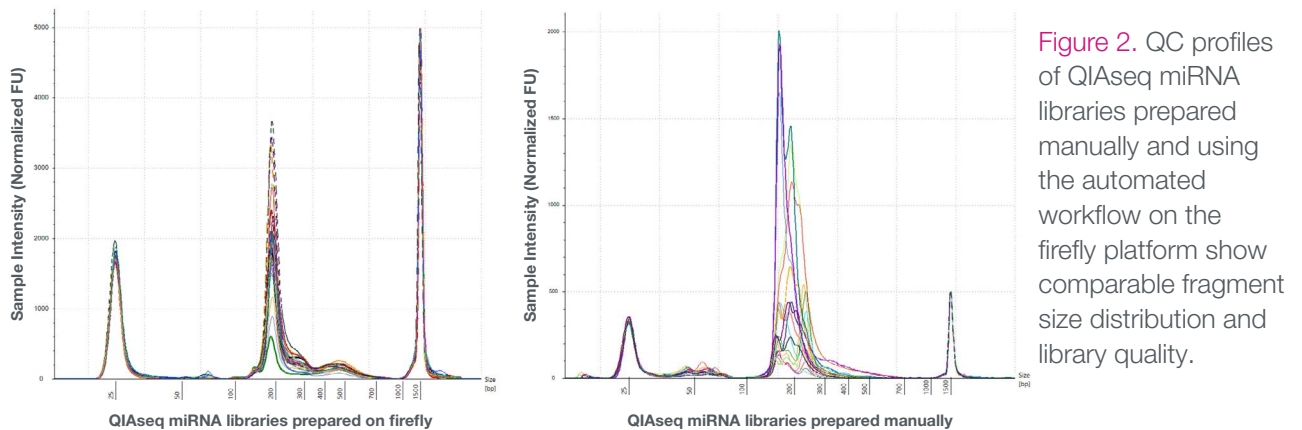
The QIAseq[®] miRNA Library Kit enables Sample to Insight, precision next-generation sequencing (NGS) of mature miRNAs on Illumina NGS instruments. This highly optimized solution facilitates both enhanced differential expression analysis using integrated Unique Molecular Indices (UMIs) and novel discovery of miRNA from cells, tissues and biofluids. Here we demonstrate the successful automation of the QIAseq miRNA workflow on the compact and intuitive [firefly[®]](#) liquid handling platform.

Workflow overview



Protocol performance

QIAseq miRNA libraries generated using automated and manual workflows*



*The data used in this application were generated by Biopolymer Facility (BPF) at Harvard Medical School.

Conclusion

firefly enables reproducible automation of the QIAseq miRNA workflow, generating libraries with consistent fragment size distribution and concentration suitable for downstream sequencing.

