



revvity

Gel-free small RNA library prep kit for Illumina[®] sequencing

NEXTFLEX[®] small RNA-seq kit v4 for Illumina[®] platforms

- Totally gel-free from miRNA or 1 ng of total RNA or more
- Completely automated protocol on the Revvity[®] Sciclone[®] G3 NGSx, Sciclone[®] G3 NGSx iQ™, and the Zephyr[®] G3 NGS workstation
- Patent pending technology optimized for challenging samples such as plasma or serum, including exosomes
- Exceptional miRNA discovery
- Ready to sequence libraries in less than 6 hours
- Multiplexing of up to 384 samples
- An automation protocol is available



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

Gel-free small RNA library prep with reduced bias

The NEXTFLEX® Small RNA-Seq Kit v4 uses patent-pending technology to provide a completely gel-free small RNA library preparation solution for Illumina® sequencing platforms. Our approach to reducing bias involves optimization of the ligation reaction and the use of dimer reduction oligonucleotides. These changes result in decreased bias in comparison to NEXTFLEX® Small RNA-Seq Kit v3 and standard protocols.

NEXTFLEX® Small RNA-Seq Kit v4 allows preparation of libraries ready for sequencing with a simplified workflow in approximately 5 hours, with as little input as 1 ng of total RNA. Libraries prepared with this kit have a higher proportion of mapping reads and unique miRNAs than leading competitors (Figure 1 and 2).

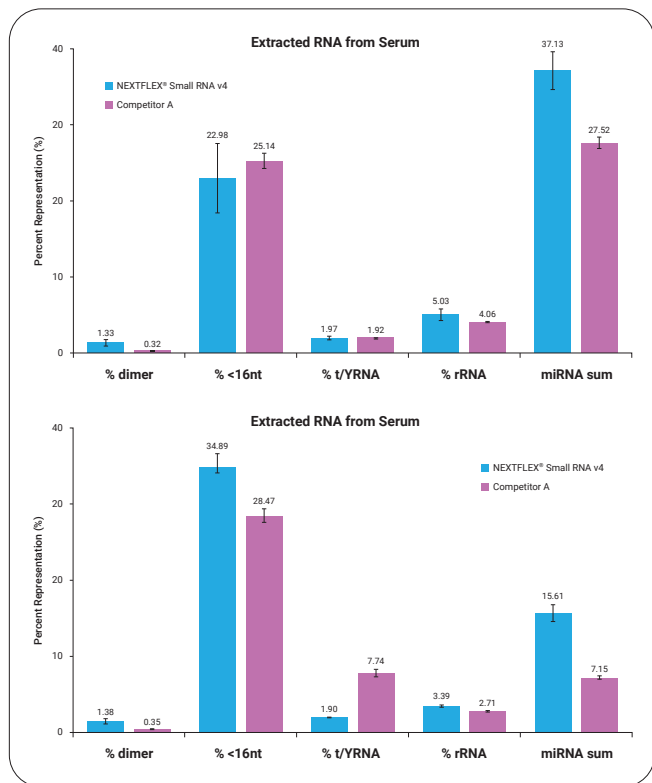


Figure 1

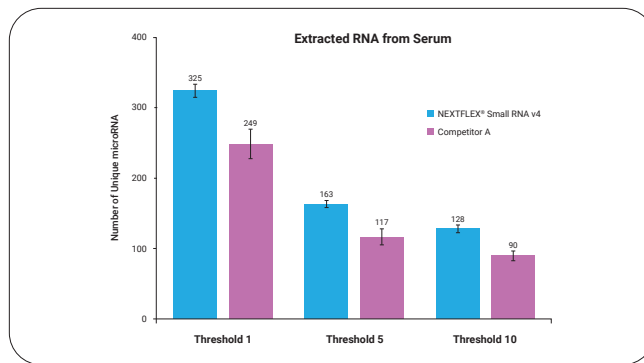


Figure 2

Maximize miRNA mapped reads with NEXTFLEX® blood miRNA blockers

The NEXTFLEX® Blood miRNA blockers deplete common miRNA in blood and plasma such as miR-486-5p, miR-92a-3p and miR-451a. These constitute 50-70% of the small RNA present in those samples and usually are not of interest for researchers.

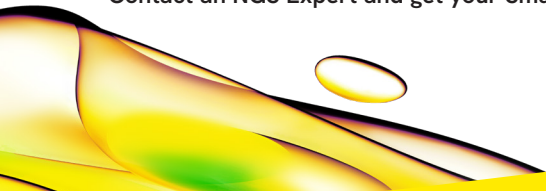
Small RNA-seq library prep optimized for automation

By eliminating all gel-based steps even for low input samples, the NEXTFLEX® Small RNA-Seq Kit V4 is optimized for easy migration onto automated liquid handling platforms. Currently methods are available for the Revvity Sciclone® G3 NGSx, Sciclone G3 NGSx iQ™, and the Zephyr® G3 NGS workstation.

Illumina® small RNA-seq multiplexing with UDI barcodes

Up to 384 UDI barcoded primers are included in each kit for multiplexing. These UDIs have been redesigned allowing libraries prepared with this kit can be run on any Illumina® sequencing platform, including NextSeq® and Novaseq® sequencers using standard Illumina® sequencing primers.

Contact an NGS Expert and get your Small RNA project started today!



revvity