



# Simplify & reduce the cost of hybridization-based target capture with universal blockers.

## NEXTFLEX® universal blockers

- Effective in blocking single-index or dual-index libraries
- Compatible with all NEXTFLEX® barcodes and all other ligation-based and tagmentation-based workflows for Illumina® sequencing
- Simplifies and reduces the cost of target capture without sacrificing performance

## Improve the number of on-target reads

During target enrichment, adapters can interact with the complementary adapter sequence strand of another library molecule. The NEXTFLEX® Universal Blockers are designed

to prevent concatemerization of library molecules. By blocking this complementary adapter sequence interaction, the number of on-target reads improves dramatically and the depth of enrichment increases.

## One set of blockers for all applications

The NEXTFLEX® Universal Blockers are designed for use with indexed adapters for Illumina® sequencing platforms. They are effective in blocking single-indexed or dual-indexed libraries, without being affected by index length or the presence of a unique molecular index (UMI).

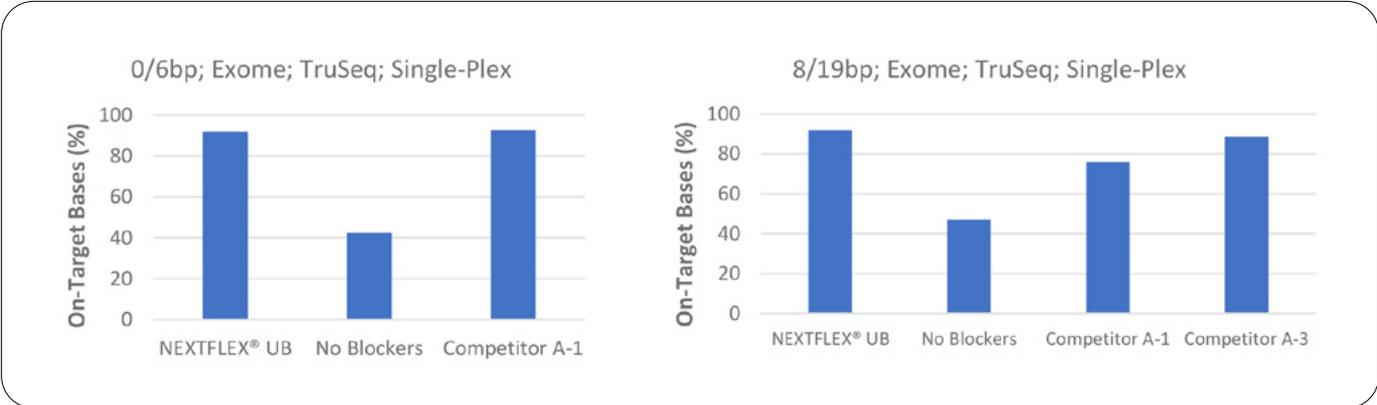


Figure1: NEXTFLEX® Universal Blockers (UB) work from 0-19 bp unblocked regions (barcodes and/or UMIs). Competitor A1- and A-3 are two separate products from Competitor A.

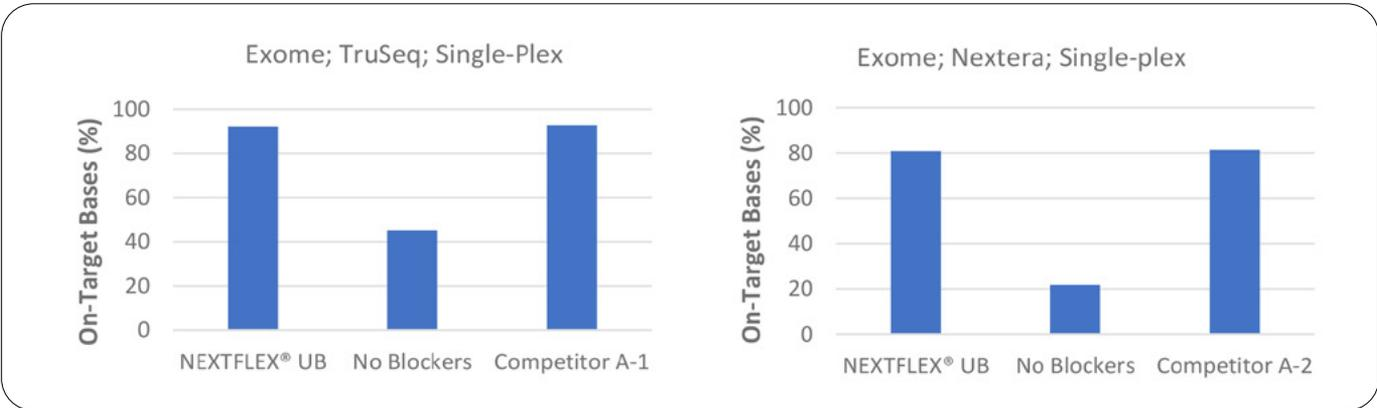


Figure2: NEXTFLEX® Universal Blockers can block Illumina® TruSeq™ or Nextera™-style libraries. Competitor A1- and A-2 are two separate products from Competitor A.

## Compatible with multiple workflows

The NEXTFLEX® Universal Blockers are compatible with ligation-based workflows, such as NEXTFLEX® or Illumina® TruSeq™ library prep kits, and with any libraries prepared with the NEXTFLEX® Unique Dual Index Barcodes.

They are also compatible with tagmentation-based workflows, such as Illumina® Nextera™ library prep kits.

Contact Us to Simplify & Reduce the Cost of Your Target Capture

