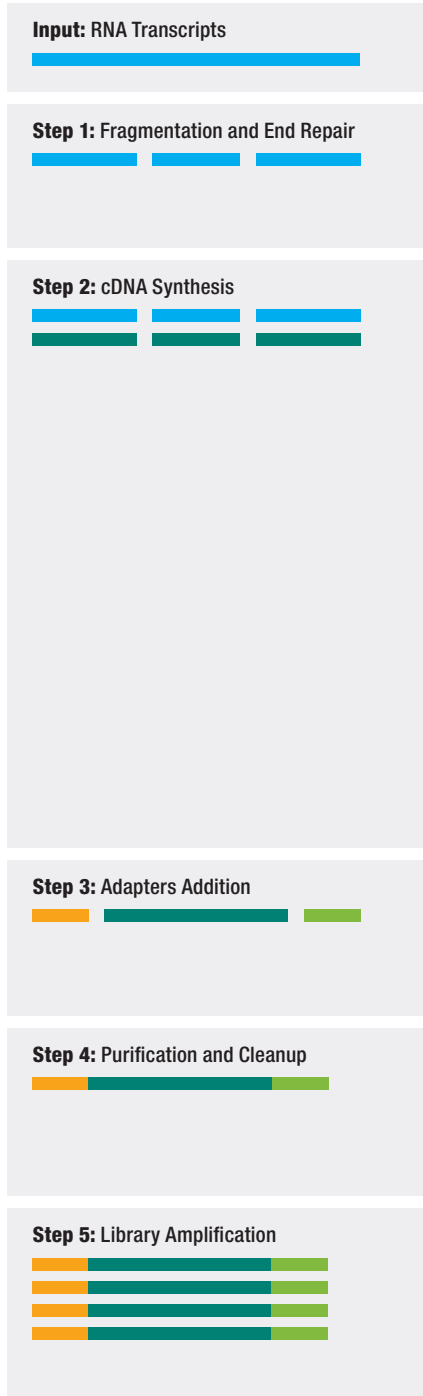


Capture More of the Transcriptome in Less Time

Traditional RNA Library Prep Workflow



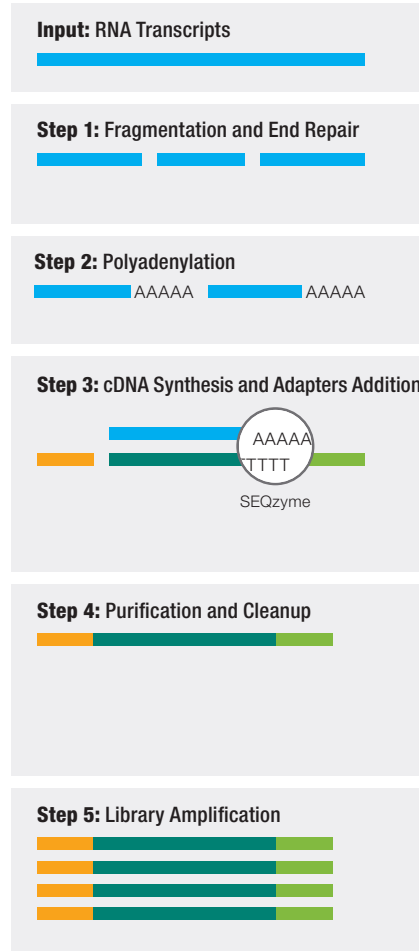
TUBES PER KIT: 15+



TOTAL TIME:

6-8
hours

SEQuoia Complete Stranded RNA Library Prep Kit Workflow



TUBES PER KIT: 7



TOTAL TIME:

<4
hours

ADVANTAGES

Use a broad range of input quantities (100 pg to 1 µg)

Avoid enzymatic bias by using a non-enzymatic fragmentation method

Capture all RNA biotypes, regardless of size or abundance

Eliminate reverse transcriptase and ligase bias using a unique continuous synthesis reaction

Maintain >99% strandedness

Increase throughput with the automation-friendly, magnetic bead-based purification protocol

Reduce duplication rate with fewer PCR cycles and better library yield

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.

Bulletin 3268 Ver A 21-0697 1021