

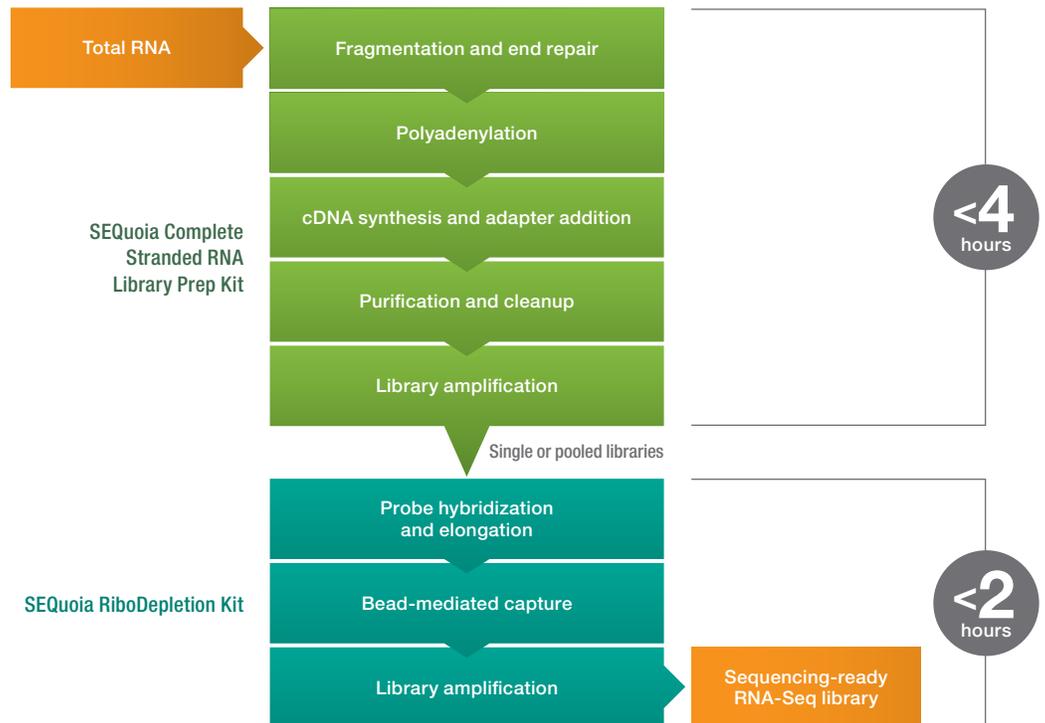
Explore the SEQuoia Workflow for Revolutionized RNA Sequencing

Profile the Whole Transcriptome with an Easy and Streamlined Workflow

- Unique patented enzymatic technology captures both long and short RNAs
- Superior performance captures more RNA fragments from low-input and low-quality samples
- Continuous synthesis reaction minimizes pipetting steps and simplifies the workflow
- Cost-effective workflow enables ribodepletion on pooled libraries

Overcome the limitations imposed by low-quantity and degraded RNA samples and capture more of the transcriptome using the SEQuoia workflow. Featuring the latest advancement in RNA sequencing (RNA-Seq) library preparation chemistry, this workflow captures and retains both long and short RNA fragments and is ideal for use with challenging samples.

The streamlined protocol can be completed in less than 6 hours.



SEQuoia Complete Stranded RNA Library Prep Kit



A high-performance RNA–Seq library prep kit that efficiently captures all RNA subtypes in a single library, even from limited and low-quality samples.

- Ideal for limited or degraded RNA
- Validated with formalin-fixed paraffin-embedded and other low-quality samples
- Streamlined protocol (less than 4 hours)
- Integrated with SeqSense NGS Data Analysis Software, a bioinformatics solution
- Compatible with all Illumina® sequencing platforms

SEQuoia Dual Indexed Primers



High-quality, unique dual indexed (UDI) primer pairs designed for use with the SEQuoia Complete Stranded RNA Library Prep Kit.

- Balanced for optimal multiplexed sequencing results
- Available in 12 UDI or 96 UDI configuration

SEQuoia RiboDepletion Kit



A stand-alone, post-library prep ribosomal RNA depletion kit that minimizes the number of library constructs mapping to ribosomal RNA while retaining more of the transcriptome.

- Efficient removal of RNA-Seq library fragments derived from ribosomal RNA
- Ideal for limited or degraded RNA samples
- Compatible with most RNA library prep kits for use on Illumina sequencing platforms
- Flexible protocol enables depletion of multiple libraries in a single reaction

Ordering Information

Catalog #	Description
17005726	SEQuoia Complete Stranded RNA Library Prep Kit , 24 reactions
17005710	SEQuoia Complete Stranded RNA Library Prep Kit , 96 reactions
12011928	SEQuoia Dual Indexed Primers Set , 12 vials of unique dual indexed primers, 96 reactions
12011930	SEQuoia Dual Indexed Primers Plate , 96-well plate of unique dual indexed primers, 96 reactions
17006487	SEQuoia RiboDepletion Kit , 24 reactions

Visit bio-rad.com/NGS for more information.

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



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

