

## Explore the SEQuoia Workflow for Revolutionized RNA Sequencing

- Patented enzymatic technology captures more of the transcriptome
- Optimized reaction generates greater yields with fewer PCR cycles
- Simplified workflow minimizes pipetting steps
- Increased cost savings by multiplexing ribodepletion reactions
- Integrated bioinformatics solution conveniently processes FASTQ files

Overcome the limitations of complicated and biased library prep methods using the SEQuoia RNA-Seq library prep workflow to capture more of the transcriptome. Featuring the latest advancements in patented enzymatic technology, SEQuoia products offer superior construction of RNA-Seq libraries from low-input and low-quality samples. The SEQuoia one-tube continuous synthesis library prep reaction minimizes pipetting steps and shortens the workflow. When used in conjunction with the SEQuoia Dual Indexed Primers and the post-library SEQuoia RiboDepletion Kit, the SEQuoia workflow can greatly reduce the total cost of library preparation without compromising data integrity or quality.

SeqSense NGS Data Analysis Software is a hosted bioinformatics pipeline uniquely optimized for mapping, aligning, counting, and normalizing reads generated from SEQuoia RNA-Seq libraries. Perform secondary analysis, determine the quality of the sequencing dataset, and generate output files ready for tertiary analysis. Analysis toolkits are also available as downloadable docker containers for deployment on local infrastructure.

### Explore More with the SEQuoia Workflow

#### Library Preparation

##### SEQuoia Complete Stranded RNA Library Prep Kit

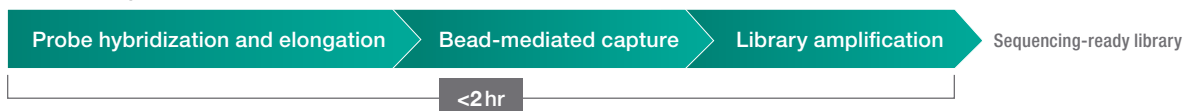


##### SEQuoia Express Stranded RNA Library Prep Kit



#### Ribodepletion

##### SEQuoia RiboDepletion Kit



**Table 1. SEQuoia Product Selection.**

	SEQuoia Complete Stranded RNA Library Prep Kit	SEQuoia Express Stranded RNA Library Prep Kit	SEQuoia Dual Indexed Primers	SEQuoia RiboDepletion Kit
Capture range	All RNA subtypes >20 bp	Long RNA fragments >200 bp	Unique dual indexed primer set optimized for performance with SEQuoia Complete and SEQuoia Express Library Prep Kits	Post-library depletion kit compatible with most RNA-Seq libraries, including those constructed using the SEQuoia Complete and SEQuoia Express Library Prep Kits
Input range	100 pg–1 µg	1 ng–1 µg		
Sample type	All sample types, including low-quality formalin-fixed paraffin-embedded (FFPE) samples	Standard good-quality sample		
Sequencing read	Single-read	Paired-end		
Protocol time	4 hr	3 hr		

### 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 FFPE and other low-quality samples
- Streamlined protocol (<4 hr)
- Integrated with SeqSense NGS Data Analysis Software, a bioinformatics solution
- Compatible with all Illumina® sequencing platforms

### SEQuoia Express Stranded RNA Library Prep Kit



A 3-tube, 3-hr library preparation solution that efficiently captures mRNA and long noncoding RNA fragments.

- Efficient and unbiased capture of mRNA and long noncoding long steps
- Integrated bioinformatics solution

### 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	<b>SEQuoia Complete Stranded RNA Library Prep Kit</b> , 24 reactions
17005710	<b>SEQuoia Complete Stranded RNA Library Prep Kit</b> , 96 reactions
12017297	<b>SEQuoia Express Stranded RNA Library Prep Kit</b> , 24 reactions
12017265	<b>SEQuoia Express Stranded RNA Library Prep Kit</b> , 96 reactions
12011928	<b>SEQuoia Dual Indexed Primers Set</b> , 12 vials of unique dual indexed primers, 96 reactions
12011930	<b>SEQuoia Dual Indexed Primers Plate</b> , 96-well plate of unique dual indexed primers, 96 reactions
17006487	<b>SEQuoia RiboDepletion Kit</b> , 24 reactions

Visit [bio-rad.com/NGS](https://www.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. © 2022 Bio-Rad Laboratories, Inc.



**Bio-Rad  
Laboratories, Inc.**

---

Life Science  
Group

**Website** [bio-rad.com](http://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

