



January 2026

QIAseq[®] FastSelect[®] Handbook

Removal of unwanted RNA, including ribosomal RNA (rRNA), globin mRNA, and mitochondrial genes for RNA-seq library preparation

Table of Contents

- Kit Contents 5
- Shipping and Storage 7
- Intended Use 8
- Safety Information 9
- Quality Control 10
- Introduction 11
 - Principle and procedure 13
 - QIAseq FastSelect: Design 14
 - QIAseq FastSelect –rRNA HMR: rRNAs removed and species covered 14
 - QIAseq FastSelect –Globin HMR: Globin genes removed and species covered 14
 - QIAseq FastSelect –rRNA Plant: rRNAs removed and species covered 15
 - QIAseq FastSelect –rRNA Yeast: rRNAs removed 15
 - QIAseq FastSelect –rRNA Worm: rRNAs removed 15
 - QIAseq FastSelect –rRNA Fish: rRNAs removed 15
 - QIAseq FastSelect –rRNA Fly: rRNAs removed 16
 - QIAseq FastSelect –Mitochondrial Genes RNAs removed 16
 - QIAseq FastSelect: Sample type and short-read RNA-seq library kit compatibility .. 16
 - QIAseq FastSelect: Sample type and long-read compatibility 17
- Important Notes 18
- Protocol: FastSelect –rRNA, –Globin mRNA, or –Mitochondrial Genes with the QIAseq Stranded Total RNA Library Kits 20
 - Important points before starting 20
 - Procedure 21
- Protocol: FastSelect –Globin with the QIAseq Stranded mRNA Select Kit 24
 - Important points before starting 24
 - Procedure 24
- Protocol: FastSelect –rRNA and/or –Globin with Illumina Stranded Total RNA Prep 27
 - Important points before starting 27
 - Procedure 27

Protocol: FastSelect –rRNA and/or –Globin with Illumina Stranded mRNA Prep	30
Important points before starting	30
Procedure	30
Protocol: FastSelect –rRNA and/or –Globin with TruSeq Stranded Library Preparation	33
Important points before starting	33
Procedure	33
Protocol: FastSelect –Globin with the TruSeq Stranded mRNA Library Preparation	36
Important points before starting	36
Procedure	36
Protocol: FastSelect –rRNA and/or –Globin with the NEBNext Ultra II Directional Library Prep Kit	38
Important points before starting	38
Procedure	38
Protocol: FastSelect –Globin with the NEBNext Poly(A) mRNA Magnetic Isolation Module and NEBNext Ultra II Directional Library Prep Kit	41
Important points before starting	41
Procedure	41
Protocol: FastSelect –rRNA and/or –Globin with the KAPA RNA HyperPrep Kit	44
Important points before starting	44
Procedure	44
Protocol: FastSelect –Globin with the KAPA mRNA HyperPrep Kit	49
Important points before starting	49
Procedure	49
Troubleshooting Guide	52
Appendix A: Species Compatibility	53
QIAseq FastSelect –rRNA HMR and –Globin species compatibility	53
QIAseq FastSelect –rRNA Plant: Types of rRNAs that are removed and species	55
QIAseq FastSelect –Mitochondrial Genes Kit	56
Appendix B: FastSelect –rRNA, –Globin mRNA, and Mitochondrial Genes with the QIAseq Stranded RNA Lib Kit	57
Important points before starting	57
Procedure	57

Ordering Information 62

Document Revision History 65

Kit Contents

Human/Mouse/Rat (HMR)

QIAseq FastSelect –rRNA HMR Kit	(24)	(96)	(384)
Catalog no.	334386	334387	334388
Number of reactions	24	96	384
QIAseq FastSelect –rRNA HMR	3 × 12 µL	120 µL	4 × 120 µL
QIAseq FastSelect –Globin Kit	(24)	(96)	(384)
Catalog no.	334376	334377	334378
Number of reactions	24	96	384
QIAseq FastSelect –Globin	3 × 12 µL	120 µL	4 × 120 µL
QIAseq FastSelect –rRNA/Globin Kit	(24)	(96)	(384)
Catalog no.	335376	335377	335378
Number of reactions	24	96	384
QIAseq FastSelect –rRNA HMR	3 × 12 µL	120 µL	4 × 120 µL
QIAseq FastSelect –Globin	3 × 12 µL	120 µL	4 × 120 µL

Mitochondrial genes

QIAseq FastSelect –Mitochondrial Genes Kit	(24)	(96)
Catalog no.	333282	333286
Number of reactions	24	96
QIAseq FastSelect –Mitochondrial Genes	36 µL	120 µL

Plant

QIAseq FastSelect –rRNA Plant Kit	(24)	(96)	(384)
Catalog no.	334315	334317	334319
Number of reactions	24	96	384
QIAseq FastSelect –rRNA Plant	3 × 12 µL	120 µL	4 × 120 µL

Yeast

QIAseq FastSelect –rRNA Yeast Kit	(24)	(96)	(384)
Catalog no.	334215	334217	334219
Number of reactions	24	96	384
QIAseq FastSelect –rRNA Yeast	3 × 12 µL	120 µL	4 × 120 µL

Worm

QIAseq FastSelect –rRNA Worm Kit	(24)	(96)
Catalog no.	333242	333245
Number of reactions	24	96
QIAseq FastSelect –rRNA Worm	3 × 12 µL	120 µL

Fish

QIAseq FastSelect –rRNA Fish Kit	(24)	(96)
Catalog no.	333252	333255
Number of reactions	24	96
QIAseq FastSelect –rRNA Fish	3 × 12 µL	120 µL

Fly

QIAseq FastSelect –rRNA Fly Kit	(24)	(96)
Catalog no.	333262	333265
Number of reactions	24	96
QIAseq FastSelect –rRNA Fly	3 × 12 µL	120 µL

Shipping and Storage

QIAseq FastSelect is shipped on blue ice or dry ice. Upon receipt, all components should immediately be stored in a constant-temperature freezer at -30°C to -15°C . Under these conditions, the components are stable, without showing any reduction in performance and quality, until the date indicated on the box label.

Intended Use

All QIAseq FastSelect products are intended for molecular biology applications. These products are not intended for the diagnosis, prevention, or treatment of a disease.

All due care and attention should be exercised in the handling of the products. We recommend all users of QIAGEN® products to adhere to the NIH guidelines that have been developed for recombinant DNA experiments, or to other applicable guidelines.

Safety Information

When working with chemicals, always wear a suitable lab coat, disposable gloves, and protective goggles. For more information, please consult the appropriate safety data sheets (SDSs). These are available online in convenient and compact PDF format at www.qiagen.com/safety, where you can find, view, and print the SDS for each QIAGEN kit and kit component.

Quality Control

In accordance with QIAGEN's ISO-certified Quality Management System, each lot of QIAseq FastSelect is tested against predetermined specifications to ensure consistent product quality.

Introduction

RNA-focused next-generation sequencing (NGS) enables a thorough investigation of both coding and noncoding RNAs. While performing RNA library preparation, significantly over-represented RNAs such as ribosomal RNA (rRNA), globin messenger RNA (mRNA) and mitochondrial genes may be avoided to facilitate optimal read allocation. QIAseq FastSelect is a breakthrough technology that rapidly and efficiently removes both cytoplasmic and mitochondrial rRNA, globin mRNA and mitochondrial genes during NGS library preparation starting with 1 ng to 1 µg of total RNAs. QIAseq FastSelect is available for multiple species and will work on non-tested related species based on sequence homology.

Cytoplasmic and mitochondrial rRNA comprise more than 80% of the total RNA in common biological samples, while globin mRNA is significantly overrepresented in whole-blood samples. As a result, commercial solutions exist to either enrich for poly(A)⁺ RNAs or deplete rRNA and/or globin. Existing strategies to remove rRNA and highly expressed mRNAs rely on probe-based depletion or enzymatic (RNase H) removal strategies. Unfortunately, these depletion methods are tedious, often taking more than 2 hours with extensive sample handling steps and are unable to work with formalin-fixed paraffin-embedded (FFPE) or fragmented samples.

Using the appropriate QIAseq FastSelect, the removal of rRNA, globin mRNA, and mitochondrial genes from a variety of species types can be accomplished in just 14 minutes.

QIAseq FastSelect works with existing RNA-seq workflows for the removal of unwanted RNAs during the reverse transcription step of NGS library preparation and is compatible with a broad range of commercially available RNA-seq library preparation kits. In a single step, QIAseq FastSelect removes up to 99% of unwanted cytoplasmic and mitochondrial rRNA, globin mRNA, and mitochondrial genes from a variety of species types (Table 1) in an assortment of RNA libraries (Table 2).

Table 1. Summary of available QIAseq FastSelect –rRNA and globin species

Specific species	rRNA	Globin	Mitochondrial genes
Human, Mouse, and Rat (HMR)	Yes	Yes	Yes
Plant	Yes	No	No
Yeast	Yes	No	No
Worm	Yes	No	No
Fish	Yes	No	No
Fly	Yes	No	No

Table 2. QIAseq FastSelect Kit compatibility*

Vendor	Kit	Cat. no.	Total RNA range
QIAGEN	QIAseq Stranded RNA Lib Kit UDI	180450, 180451, 180452, 180453, 180454	100 ng to 1 µg
	QIAseq Stranded mRNA Lib Kit UDI	180440, 180441, 180442, 180443, 180445	100 ng to 1 µg
	QIAseq Low Input RNA Library Kits	334202, 334205, 334207, 334208	1 ng to 1 µg
	QIAseq FastSelect RNA Library Kits	334262, 334222, 334242, 334272, 334302, 334232, 334252, 334292, 334282	1 ng to 1 µg
Illumina®	Illumina Stranded Total RNA Prep	20040525, 20040529	1 ng to 1 µg
	Illumina Stranded mRNA Prep	20040532, 20040534	25 ng to 1 µg total RNA
	TruSeq® Stranded Lib Prep	20020596, 20020597	100 ng to 1 µg
	TruSeq Stranded mRNA Lib Prep	20020594, 20020595	100 ng to 1 µg total RNA
	TruSeq RNA Library prep Kit v2	RS-122-2001, RS-122-2002	10–400 ng mRNA 100 ng to 1 µg total RNA

Table 2. QIAseq FastSelect Kit compatibility* (continued)

Vendor	Kit	Cat. no.	Total RNA range
New England Biolabs	NEBNext® Ultra™ II Directional	E7760S, E7760L	5 ng to 1 µg
Kapa® Biosystems	KAPA RNA HyperPrep	KK8540, KK8541	25 ng to 1 µg
	KAPA mRNA HyperPrep	KK8580, KK8581	50 ng to 1 µg intact total RNA
MGI Tech Co	MGIEasy Fast RNA Library Prep Set	940-000890-00, 940-000889-00, 940-001835-00, 940-000961-00	10 ng to 1 µg
Watchmaker Genomics	Watchmaker RNA Library Prep Kit	7K0078-024, 7K0078-096, 7K0078-384	1 ng to 100 ng
	Watchmaker mRNA Library Prep Kit	7BK0001-024, 7BK0001-096, 7BK0001-384	2.5 ng to 1000 ng
Takara	SMART-Seq Total RNA Mid Input	635049, 635050, 635048	1 ng to 100 ng

* QIAseq FastSelect is compatible with almost any stranded RNA library prep kit. Ensure to incorporate the depletion incubation steps prior reverse transcription. For questions regarding kits that are not listed, please contact QIAGEN Technical Services.

Principle and procedure

QIAseq FastSelect is designed for fast and selective removal of cytoplasmic and mitochondrial rRNA, globin mRNA, and mitochondrial genes from total RNA during NGS RNA library preparation. The simple one-step protocol calls for the appropriate FastSelect reagent to be directly combined with total RNA (1 ng to 1 µg) and any reaction buffers necessary for Mg²⁺ heat-dependent RNA fragmentation. After the optional heat fragmentation is performed, the reaction is gradually cooled to room temperature (15–25°C). The remaining library preparation–specific steps are then followed without the need for any additional enzymatic cleanups or RNA depletion steps. QIAseq FastSelect has been designed to work equally well with high-quality RNA or highly fragmented samples.

QIAseq FastSelect: Design

QIAseq FastSelect –rRNA HMR: rRNAs removed and species covered

QIAseq FastSelect –rRNA HMR has been designed to remove cytoplasmic (5S, 5.8S, 18S, and 28S), mitochondrial (12S and 16S), rRNA from human, mouse, and rat. In addition to human, mouse, and rat samples, QIAseq FastSelect –rRNA HMR has been tested with a variety of other RNA samples (Appendix A: Species Compatibility). Depending on rRNA sequence homology, QIAseq FastSelect –rRNA HMR will work on additional species.

QIAseq FastSelect –Globin HMR: Globin genes removed and species covered

QIAseq FastSelect –Globin HMR has been designed to remove all subunits of the hemoglobin family in human, mouse, and rat. In human specifically: HBA1, HBA2, HBB, and HBD; fetal: HBG1, HBG2, and HBQ1; embryonic: HBE1 and HBZ. In mouse specifically: Hba-a1, Hba-a2, Hbb-b1, Hbb-b2, Hbb-bs, and Hbb-bt; embryonic: Hbb-y, Hbb-bh1, and Hba-x. In rat specifically: Hba1, Hba-a1, Hba2, Hbb-b1, Hbb, Hbg1, Hbe1, Hbe2, Hbz, and Hbq1b. In addition to human, mouse, and rat samples, QIAseq FastSelect –Globin may work on species beyond human, mouse, and rat depending on globin sequence homology.

Important: When working with total RNA blood samples, QIAseq FastSelect –rRNA HMR must be used in combination with QIAseq FastSelect –Globin. When working with mRNA-enriched samples, only QIAseq FastSelect –Globin needs to be used.

QIAseq FastSelect –rRNA Plant: rRNAs removed and species covered

QIAseq FastSelect –rRNA Plant has been designed to remove cytoplasmic (5.8S, 18S, and 25S), mitochondrial (5S, 18S, and 26S), and chloroplast (4.5S, 5S, 16S, and 23S) rRNA from *Arabidopsis thaliana* and *Arabidopsis lyrata*. In addition to *Arabidopsis* total RNA, QIAseq FastSelect –rRNA Plant has been tested with a variety of other RNA samples (Appendix A: Species Compatibility, page 53). Depending on rRNA sequence homology, QIAseq FastSelect –rRNA Plant will work on additional plant species.

QIAseq FastSelect –rRNA Yeast: rRNAs removed

QIAseq FastSelect –rRNA Yeast has been designed to remove cytoplasmic (35S made up of 25S 18S & 5.8S, 25S, 18S, 5.8S, and 5S) and mitochondrial (21S, 15S, ACI60_gr01 [large subunit], and ACI60_gr02 [small subunit]) rRNA based on sequence information from *Saccharomyces cerevisiae* and *Schizosaccharomyces pombe*. Depending on rRNA sequence homology, QIAseq FastSelect –rRNA Yeast will work on additional yeast species.

QIAseq FastSelect –rRNA Worm: rRNAs removed

QIAseq FastSelect –rRNA Worm has been designed to remove 5S, 5.8S, 18S, 26S, and mitochondrial rRNA from *Caenorhabditis elegans* (roundworm). Depending on rRNA sequence homology, QIAseq FastSelect –rRNA Worm will work on additional worm species.

QIAseq FastSelect –rRNA Fish: rRNAs removed

QIAseq FastSelect –rRNA Fish has been designed to remove all annotated rRNAs in RefSeq and Ensembl for *Danio rerio* (Zebrafish) and *Salmo salar* (Atlantic salmon). Depending on rRNA sequence homology, QIAseq FastSelect –rRNA Fish will work on additional fish species.

QIAseq FastSelect –rRNA Fly: rRNAs removed

QIAseq FastSelect –rRNA Fly has been designed to remove rRNAs in 2S, 5S, 5.8S, 18S, 28S, and mitochondrial rRNA based on sequence information from *Drosophila melanogaster*. Depending on rRNA sequence homology, QIAseq FastSelect –rRNA Fly will work on additional fly species.

QIAseq FastSelect –Mitochondrial Genes RNAs removed

QIAseq FastSelect –Mitochondrial Genes has been designed to remove human, mouse, and rat mitochondrial genes:

- **Human (hg-38)** — MT-RNR1, MT-RNR2, MT-ND1, MT-ND2, MT-CO1, MT-CO2, MT-ATP8, MT-ATP6, MT-CO3, MT-ND3, MT-ND4L, MT-ND4, MT-ND5, MT-ND6, and MT-CYB
- **Mouse (mm10)** — mt-Rnr1, mt-Rnr2, mt-Nd1, mt-Nd2, mt-Co1, mt-Co2, mt-Atp8, mt-Atp6, mt-Co3, mt-Nd3, mt-Nd4l, mt-Nd4, mt-Nd5, mt-Nd6, and mt-Cytb
- **Rat (rn6)** — AY172581.9, AY172581.24, Mt-nd1, Mt-nd2, Mt-co1, Mt-co2, Mt-atp8, Mt-atp6, Mt-co3, Mt-nd3, Mt-nd4l, Mt-nd4, Mt-nd5, Mt-nd6, and Mt-cyb

QIAseq FastSelect: Sample type and short-read RNA-seq library kit compatibility

QIAseq FastSelect has been designed to be compatible with total RNA- and mRNA-enriched samples isolated from cells, fresh/frozen tissue, FFPE tissue, whole blood, and serum/plasma samples, including exosomes. The QIAseq FastSelect reagent and protocol has been tested with a variety of commercially available stranded RNA library preparation kits from QIAGEN, Illumina, NEB®, Takara Bio®, Watchmaker®, and Roche®. For questions regarding specific protocols for kits that are not detailed in the handbook, please contact QIAGEN Technical Services.

QIAseq FastSelect: Sample type and long-read compatibility

QIAseq FastSelect has been successfully used with high quality RNA for cDNA sequencing on the Oxford Nanopore. For recommendations on long-read sequencing, please contact QIAGEN Technical Services.

As a general recommendation, you can use FastSelect depletion before the RT step. To maintain fragment length, reduce the incubation time starting from the step at 65–70°C. This will allow efficient depletion without increasing RNA fragmentation.

Important Notes

- QIAGEN developed a specific low-input RNA library kit for use with FastSelect when starting with 1 ng or more of total RNA. Please see the QIAseq FastSelect RNA library kit for use with this product line (www.qiagen.com/HB-3152).
- For QIAseq FastSelect –5s/16s/23s, please see *QIAseq FastSelect –5S/16S/23S Handbook* (www.qiagen.com/HB-2695).
- We highly recommend DNase treatment (on-column or in-solution) of total RNA samples before starting your RNA-seq library preparation.
- It is not possible to test the efficiency of the FastSelect reaction by running a portion of the eluate from the bead cleanup on a TapeStation®, Fragment Analyzer™, etc. FastSelect works by inhibiting reverse transcription of rRNA, which does not occur until the first-strand synthesis reaction during library prep.
- QIAseq FastSelect is an inline solution for the removal of unwanted rRNAs during NGS library preparation. The total RNA input is defined by the range of the RNA library kit used. For example, the QIAseq Stranded Total RNA Lib Kit (cat. no. 180743 or 180745) has a total RNA input range of 100 ng to 1 µg. As a result, you would start with 100 ng to 1 µg into the FastSelect reaction. For lower input samples, the QIAseq Low Input RNA Library Kit (cat no. 334202, 334205, 334207, or 334208) can accommodate samples starting at 1 ng of total RNA.
- QIAseq FastSelect products are all compatible with each other and can be combined to remove additional RNAs. For example, when working with mixed samples, QIAseq FastSelect products are fully compatible with QIAseq FastSelect –5S/16S/23S (which is used to remove bacterial rRNA). For relevant protocols, please refer to the *QIAseq FastSelect –5S/16S/23S Handbook* (www.qiagen.com/HB-2695) or the *QIAseq FastSelect Custom Handbook* (www.qiagen.com/HB-2917).

- The rRNA removal imparted by QIAseq FastSelect is extremely robust, especially when compared to other methods. We recommend preparing libraries and using the standard protocol for library preparation unless specifically noted in the handbook.
- If the yield of the library is less than with other RNA removal methods, this is often caused by the increased removal of RNA imparted by the QIAseq FastSelect method and is normal, especially when combining multiple QIAseq FastSelect depletion reagent in one library preparation. In our experience, adding two cycles of library amplification is usually sufficient to increase library yield for all downstream quantification and sequencing applications.
- Depending on the RNA-seq kit and RNA input amounts, adapter–dimers may be observed. If this happens, we recommend performing a second bead-based cleanup reaction of the final library.
- For additional RNA library kit compatibility and protocol support, please contact QIAGEN Technical Services (www.qiagen.com/support/technical-support).

Protocol: FastSelect –rRNA, –Globin mRNA, or –Mitochondrial Genes with the QIAseq Stranded Total RNA Library Kits

Important points before starting

- This protocol is for combining two different FastSelect panels with the QIAseq Stranded RNA Library Kit in a single sample. For combining three or more panels, please see Appendix B: FastSelect –rRNA, –Globin mRNA, and Mitochondrial Genes with the QIAseq Stranded RNA Lib Kit.
- The QIAseq Stranded RNA Lib Kit (cat. no. 180450, 180451, 180452, 180453, or 180454) is required for use with this protocol.
- This protocol has been tested with 100 ng to 1 µg of total RNA.
- When removing globin, two additional cycles of Hi-Fi library amplification need to be performed. Similarly, when removing rRNA and mitochondrial genes, two additional cycles of Hi-Fi library amplification may be required to adjust the losses of library yields due to the removal of a high number of transcripts.
- Refer to the *QIAseq Stranded RNA Library Kit Handbook* available at www.qiagen.com/HB-2465

Procedure

1. Thaw total RNA on ice. Gently mix, briefly centrifuge to collect residual liquid from the sides of the tubes, and return to ice.
2. Prepare the reagents required for the RNA fragmentation and the appropriate QIAseq FastSelect rRNA, globin, and mitochondrial gene removal.
 - a. Thaw 5× RT Buffer, nuclease-free water from the QIAseq stranded kit, and the tube(s) from the appropriate QIAseq FastSelect kit at room temperature.
 - b. Mix by vortexing and then briefly centrifuge.
3. On ice, prepare the fragmentation/RNA Removal reaction according to Table 3. Briefly centrifuge, mix by pipetting up and down 10 times, and centrifuge briefly again.

Note: If setting up more than one reaction, prepare a volume of master mix that is 10% greater than what is required for the total number of reactions.

If omitting FastSelect, replace with water to keep the final reaction volumes consistent.

Table 3. Setup of fragmentation/RNA removal reactions

Component	Volume/reaction
Total RNA (100 ng to 1 µg)	Variable
RT Buffer, 5x*	8 µL
QIAseq FastSelect –rRNA†	1 µL
QIAseq FastSelect –Globin or QIAseq FastSelect –Mitochondrial Genes†	1 µL
ERCC Control‡	Optional
Nuclease-free water	Variable
Total volume	37 µL

* From QIAseq Stranded Total RNA Lib Kit.

† Choose the appropriate QIAseq FastSelect based on the sample type and application. Replace FastSelect with water to keep the final volumes consistent. If combining more than two panels together, see Appendix B: FastSelect –rRNA, –Globin mRNA, and Mitochondrial Genes with the QIAseq Stranded RNA Lib Kit.

‡ ERCC Control RNA can be added according to the concentrations specified by the manufacturer. If added, the total fragmentation/RNA Removal reaction volume should remain 37 µL.

4. Incubate as described in Table 4, according to input RNA quality and desired insert size.

Table 4. Combined QIAseq fragmentation and FastSelect hybridization protocol

Input RNA quality	Step	Insert size approx. 150–250 bp	Insert size approx. 350 bp
High quality (RIN >9)	1 *	15 min at 95°C	3 min at 95°C
Moderate quality (RIN 5–6)	1 *	10 min at 95°C	3 min at 95°C
FFPE or degraded sample (RIN <3)	1 *	No fragmentation†	No fragmentation†
Steps 2–9 are performed regardless of input RNA quality. They need to be performed whether the RNA is high quality, moderate quality, FFPE, or degraded.	2	2 min at 75°C	2 min at 75°C
	3	2 min at 70°C	2 min at 70°C
	4	2 min at 65°C	2 min at 65°C
	5	2 min at 60°C	2 min at 60°C
	6	2 min at 55°C	2 min at 55°C
	7	2 min at 37°C	2 min at 37°C
	8	2 min at 25°C	2 min at 25°C
	9	Hold at 4°C	Hold at 4°C

* Choose one option for the time on step 1 according to the input RNA quality and desired insert size.

† Also suitable for exosomal RNA or RNA of other origin with a size between 80 and 500 bp.

Important: Regardless of time and temperature chosen in step 1, steps 2–9 must be performed.

5. Refer to the *QIAseq Stranded RNA Library Kit Handbook* and immediately proceed to “Protocol: First-strand Synthesis”.
6. Follow the *QIAseq Stranded RNA Library Kit Handbook* to perform all remaining library construction steps.

IMPORTANT When removing globin, two additional cycles of Hi-Fi library amplification need to be performed.



Protocol: FastSelect –Globin with the QIAseq Stranded mRNA Select Kit

Important points before starting

- The QIAseq Stranded mRNA Select Kit (cat. no. 180773 or 180775) is required for use with this protocol.
- **Important:** When removing globin, two additional cycles of Hi-Fi Library Amplification need to be performed.
- Refer to the *QIAseq Stranded mRNA Library Kit Handbook* available at www.qiagen.com/HB-2464

Procedure

1. From the *QIAseq Stranded mRNA Library Kit Handbook*, perform “Protocol: mRNA Enrichment” with the recommended amount of total RNA input (100 ng to 1 µg). Ultimately elute the enriched mRNA in 10 µL. When starting the mRNA enrichment with higher input (e.g., 500 ng) and when required to run replicate or multiple analysis with the purified mRNA, you can increase the elution volume according to the input and up to 27 µL.
2. Prepare the reagents required for the RNA fragmentation and QIAseq FastSelect –Globin removal.
 - a. Thaw 5× RT Buffer, nuclease-free water from the QIAseq stranded kit, and the QIAseq FastSelect –Globin tube from the QIAseq FastSelect kit at room temperature.
 - b. Mix by vortexing and then briefly centrifuge.

3. On ice, prepare the fragmentation/RNA Removal reaction according to Table 5. Briefly centrifuge, mix by pipetting up and down 10 times, and centrifuge briefly again.

Note: If setting up more than one reaction, prepare a volume of master mix that is 10% greater than what is required for the total number of reactions.

Table 5. Setup of fragmentation/RNA removal reactions

Component	Volume/reaction
mRNA enrichment reaction (already in tube)	27 µL
RT Buffer, 5×*	8 µL
QIAseq FastSelect –Globin	1 µL
ERCC Control†	Optional
Nuclease-free water	1 µL
Total volume	37 µL

* From QIAseq Stranded Total RNA Lib Kit.

† ERCC Control RNA can be added according to the concentrations specified by the manufacturer. If added, replace the nuclease-free water (1 µL) with ERCC.

4. Incubate as described in Table 6, according to your input RNA quality and desired insert size.

Table 6. Combined QIAseq fragmentation and FastSelect hybridization protocol

Input RNA quality	Step	Insert size approx. 150–250 bp	Insert size approx. 350 bp
High quality (RIN >9)	1 *	15 min at 95°C	3 min at 95°C
Moderate quality (RIN 5–6)	1 *	10 min at 95°C	3 min at 95°C
FFPE or degraded sample (RIN <3)	1 *	No fragmentation†	No fragmentation†
Steps 2–9 are performed regardless of input RNA quality. They need to be performed whether the RNA is high quality, moderate quality, FFPE, or degraded.	2	2 min at 75°C	2 min at 75°C
	3	2 min at 70°C	2 min at 70°C
	4	2 min at 65°C	2 min at 65°C
	5	2 min at 60°C	2 min at 60°C
	6	2 min at 55°C	2 min at 55°C
	7	2 min at 37°C	2 min at 37°C
	8	2 min at 25°C	2 min at 25°C
	9	Hold at 4°C	Hold at 4°C

* Choose one option for the time on step 1 according to the input RNA quality and desired insert size.

† Also suitable for exosomal RNA or RNA of other origin with a size between 80 and 500 bp.

Important: Regardless of time and temperature chosen in step 1, steps 2–9 must be performed.

5. Refer to the *QIAseq Stranded mRNA Library Kit Handbook* and immediately proceed to “Protocol: First-strand Synthesis.”
6. Follow the *QIAseq Stranded mRNA Library Kit Handbook* to perform all remaining library construction steps.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –rRNA and/or –Globin with Illumina Stranded Total RNA Prep

Important points before starting

- The Illumina Stranded Total RNA Prep (Illumina cat. no. 20040525 and 20040529) is required for use with this protocol.

Note: Follow the steps outlined below before proceeding to the designated step in the *Illumina Stranded Total RNA Prep Ligation Reference Guide* (1000000124514 v02). By doing this, a stranded total RNA library prep will be performed.

- **Important:** It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.

Procedure

1. Thaw the tube(s) from the appropriate QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
2. To prepare 1 ng – 1 µg of total RNA, which is required to be in a maximum volume of 6.5 µL, add QIAseq FastSelect as follows:
 - **Option 1 (remove rRNA):** Add 1 µL of the appropriate QIAseq FastSelect –rRNA
 - **Option 2 (remove globin):** Add 1 µL of QIAseq FastSelect –Globin
 - **Option 3 (remove rRNA and globin):** Add 1 µL of QIAseq FastSelect –rRNA HMR and 1 µL of QIAseq FastSelect –Globin

3. From the Illumina Stranded Total RNA Library Prep, add ELB (x μL) to bring the volume of the reaction to 8.5 μL .
4. From the Illumina Stranded Total RNA Library Prep, add 8.5 μL EPH3 to bring the volume of the reaction to 17 μL .
5. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes.
6. Incubate in a thermal cycler with a heated lid as described in Table 7.

Table 7. Combined Illumina fragmentation and FastSelect hybridization protocol

Step	Time and temperature
1	2 min at 94°C
2	2 min at 75°C
3	2 min at 70°C
4	2 min at 65°C
5	2 min at 60°C
6	2 min at 55°C
7	2 min at 37°C
8	2 min at 25°C
9	Hold at 4°C

7. Using the entire 17 μL fragmented/hybridized RNA reaction, refer to the *Illumina Stranded Total RNA Prep Ligation Reference Guide* and immediately proceed to “Synthesize First Strand cDNA.”
8. Follow the *Illumina Stranded Total RNA Prep Ligation Reference Guide* to perform all remaining library construction steps.

IMPORTANT

It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.

IMPORTANT

When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –rRNA and/or –Globin with Illumina Stranded mRNA Prep

Important points before starting

- The Illumina Stranded mRNA Prep (Illumina cat. no. 20040532 and 20040534) is required for use with this protocol.

Note: With this protocol, do not perform mRNA purification. Instead, follow the steps outlined below before proceeding to the designated step in the *Illumina Stranded mRNA Prep Ligation Reference Guide* (1000000124518 v02). By doing this, a stranded total RNA library prep will be performed.

- **Important:** It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.

Procedure

1. Thaw the tube(s) from the appropriate QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
2. To prepare 25 ng to 1 µg of total RNA, which is required to be in a maximum volume of 6.5 µL, add QIAseq FastSelect as follows:
 - **Option 1 (remove rRNA):** Add 1 µL of the appropriate QIAseq FastSelect –rRNA
 - **Option 2 (remove globin):** Add 1 µL of QIAseq FastSelect –Globin
 - **Option 3 (remove rRNA and globin):** Add 1 µL of QIAseq FastSelect –rRNA HMR and 1 µL of QIAseq FastSelect –Globin

3. From the Illumina Stranded mRNA Prep, add ELB (x μ L) to bring the volume of the reaction to 8.5 μ L.
4. From the Illumina Stranded mRNA Prep, add 8.5 μ L EPH3 to bring the volume of the reaction to 17 μ L.
5. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes.
6. Incubate in a thermal cycler with a heated lid as described in Table 8.

Table 8. Combined Illumina fragmentation and FastSelect hybridization protocol

Step	Time and temperature
1	2 min at 94°C
2	2 min at 75°C
3	2 min at 70°C
4	2 min at 65°C
5	2 min at 60°C
6	2 min at 55°C
7	2 min at 37°C
8	2 min at 25°C
9	Hold at 4°C

7. Using the entire 17 μ L fragmented/hybridized RNA reaction, refer to the *Illumina Stranded mRNA Prep Ligation Reference Guide* and immediately proceed to “Synthesize First Strand cDNA.”
8. Follow the *Illumina Stranded mRNA Prep Ligation Reference Guide* to perform all remaining library construction steps.

IMPORTANT

It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.

IMPORTANT

When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –rRNA and/or –Globin with TruSeq Stranded Library Preparation

Important points before starting

- The TruSeq Stranded mRNA Library Prep (Illumina cat. no. 20020594 or 20020595) is required for use with this protocol.

Note: With this protocol, do not perform mRNA purification. Instead, follow the steps outlined below before proceeding to “Synthesize First Strand cDNA” in the *TruSeq Stranded mRNA Reference Guide*. By doing this, a stranded, total RNA library preparation will be performed.

- This protocol has been tested with 100 ng to 1 µg of total RNA.
- **Important:** It is highly recommended to dilute the Illumina adapters two-fold compared to what is suggested in the default Illumina protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *TruSeq Stranded mRNA Reference Guide* (1000000040498).

Procedure

1. Thaw the tube(s) from the appropriate QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
2. To prepare 100 ng to 1 µg of total RNA, which is required to be in a maximum volume of 5 µL, add QIAseq FastSelect as follows:
 - **Option 1 (remove rRNA):** Add 1 µL of the appropriate QIAseq FastSelect –rRNA
 - **Option 2 (remove globin):** Add 1 µL of QIAseq FastSelect –Globin

- **Option 3 (remove rRNA and globin):** Add 1 μL of QIAseq FastSelect –rRNA HMR and 1 μL of QIAseq FastSelect –Globin

3. From the TruSeq Stranded mRNA Library Prep, add 14.5 μL FPF (when using option 1 or 2 in the previous step), or add 13.5 μL FPF (when using option 3 in the previous step), to bring the volume of the reaction to 20.5 μL .
4. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes.
5. Incubate in a thermal cycler with a heated lid as described in Table 9.

Important: Table 10 can be consulted to adjust RNA insert size. Irrespective of time at 94°C, steps 2–9 listed in Table 9 must be performed.

Table 9. Combined TruSeq Stranded fragmentation and FastSelect hybridization protocol

Step	Time and temperature
1 *	8 min at 94°C*
2	2 min at 75°C
3	2 min at 70°C
4	2 min at 65°C
5	2 min at 60°C
6	2 min at 55°C
7	2 min at 37°C
8	2 min at 25°C
9	Hold at 4°C

* The initial step at 94°C can be modified to permit longer RNA insert sizes. Refer to Table 10 for recommendations.

Note: The remaining steps 2–9 are performed regardless of the time at 94°C.

Table 10. Fragmentation time at 94°C for alternative RNA insert sizes

Time at 94°C* (min)	Range of insert length (bp)	Median insert length (bp)	Average final library size (Bioanalyzer bp)
0	130–350	200	467
1	130–310	190	439
2	130–290	185	410
3	125–250	165	366
4	120–225	160	326
8	120–210	155	309
12	115–180	140	272

* The remaining steps 2–9 from Table 9 must be performed regardless of the time at 94°C.

6. Use 17 µL of the fragmented/hybridized RNA, refer to the *TruSeq Stranded mRNA Reference Guide* and immediately proceed to “Synthesize First Strand cDNA.”

Note: From the *TruSeq Stranded mRNA Reference Guide*, the procedural step “Place the RBP plate on the magnetic stand and wait until the liquid is clear (approx. 5 min)” is not applicable.

7. Follow the *TruSeq Stranded mRNA Reference Guide* to perform all remaining library construction steps.

IMPORTANT



It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –Globin with the TruSeq Stranded mRNA Library Preparation

Important points before starting

- The TruSeq Stranded mRNA Library Prep (Illumina cat. no. 20020594 or 20020595) is required for use with this protocol.
- **Important:** It is highly recommended to dilute the Illumina adapters two-fold compared to what is suggested in the Illumina protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *TruSeq Stranded mRNA Reference Guide* (1000000040498).

Procedure

1. Using the *TruSeq Stranded mRNA Reference Guide*, purify mRNA as described under “Purify mRNA” steps 1–19 (pages 11–12 in 1000000040498).
2. Using the *TruSeq Stranded mRNA Reference Guide*, fragment mRNA as described under “Fragment mRNA” steps 1–15 (page 12 in 1000000040498).
3. Using the *TruSeq Stranded mRNA Reference Guide*, perform steps 1 and 2 under the “Procedure” section (page 13 in 1000000040498) of “Synthesize First Strand cDNA”.
4. Thaw the QIAseq FastSelect –Globin tube from the QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
5. To the 17 μ L supernatant in the CDP plate, add 1 μ L of QIAseq FastSelect –Globin.
6. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes.

7. Incubate in a thermal cycler with a heated lid as described in Table 11.

Table 11. FastSelect hybridization protocol

Step	Time and temperature
1	2 min at 75°C
2	2 min at 70°C
3	2 min at 65°C
4	2 min at 60°C
5	2 min at 55°C
6	2 min at 37°C
7	2 min at 25°C
8	Hold at 4°C

8. Refer to the *TruSeq Stranded mRNA Reference Guide* and immediately proceed to and perform step 3 under the “Procedure” section (page 13) of “Synthesize First Strand cDNA”.
9. Follow the *TruSeq Stranded mRNA Reference Guide* to perform all remaining library construction steps.

IMPORTANT



It is recommended to test dilution of the Illumina anchors two-fold compared to what is suggested in the default Illumina protocol.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –rRNA and/or –Globin with the NEBNext Ultra II Directional Library Prep Kit

Important points before starting

- The NEBNext Ultra II Directional RNA Library Prep Kit for Illumina (NEB cat. no. E7760S or E7760L) is required for use with this protocol.
- This protocol has been tested with 5 ng to 1 µg of total RNA.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* (Version 3.1).

Procedure

1. Thaw the tube(s) from the appropriate QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
2. Referring to Section 4 from the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual*, perform the following in place of steps 4.1.1 through 4.1.4:
 - a. Assemble the fragmentation and priming reaction described in Table 12 on ice in a nuclease-free tube.

Table 12. NEBNext Ultra II Stranded fragmentation and priming mix

Component	Volume/reaction (μL)
Total RNA (5 ng–1 μg)	4
(lilac) NEBNext First Strand Synthesis Reaction Buffer*	4
(lilac) Random Primers*	1
Total volume	9

* From NEBNext Ultra II Directional Library Prep Kit.

- b. To the assembled fragmentation and priming mix, add QIAseq FastSelect as follows:
 - **Option 1 (remove rRNA):** Add 1 μL of the appropriate QIAseq FastSelect –rRNA
 - **Option 2 (remove globin):** Add 1 μL of QIAseq FastSelect –Globin
 - **Option 3 (remove rRNA and globin):** Add 1 μL of QIAseq FastSelect –rRNA HMR and 1 μL of QIAseq FastSelect –Globin
- c. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes
- d. Incubate in a thermal cycler with a heated lid as described in Table 13, according to your input RNA quality.

Important: Regardless of time and temperature chosen in step 1, steps 2–9 must be performed.

Table 13. Combined NEBNext Ultra II fragmentation and FastSelect hybridization protocol

Step	Intact RNA (RIN >7)	Partially degraded RNA (RIN 2–6)
1	15 min at 94°C	7–8 min at 94°C
2	2 min at 75°C	2 min at 75°C
3	2 min at 70°C	2 min at 70°C
4	2 min at 65°C	2 min at 65°C
5	2 min at 60°C	2 min at 60°C
6	2 min at 55°C	2 min at 55°C
7	2 min at 37°C	2 min at 37°C
8	2 min at 25°C	2 min at 25°C
9	Hold at 4°C	Hold at 4°C

3. Refer to the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* and immediately proceed to “First Strand cDNA Synthesis Reaction”.

Note: “First Strand cDNA Synthesis Reaction” is chapter 4.2 in version 3.1 of the instruction manual.

4. Follow the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* to perform all remaining library construction steps.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –Globin with the NEBNext Poly(A) mRNA Magnetic Isolation Module and NEBNext Ultra II Directional Library Prep Kit

Important points before starting

- The NEBNext Poly(A) mRNA Magnetic Isolation Module (NEB cat. no. E7490) is required for use with this protocol.
- The NEBNext Ultra II Directional RNA Library Prep Kit for Illumina (NEB cat. no. E7760S or E7760L) is required for use with this protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* (Version 3.1).

Procedure

1. Thaw the FastSelect –Globin tube from the QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
2. Referring to section 1 from the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual*, perform section 1.1 and section 1.2, steps 1.2.1 through 1.2.36 as indicated.
3. In place of steps 1.2.37 through 1.2.40, perform the following:
 - a. Incubate the sample in a thermal cycler (with the heated lid set at 105°C) for 15 min at 94°C but do not cool the sample to 4°C.

- b. Immediately after the 94°C fragmentation has been completed, quickly spin down the tube in a microcentrifuge to collect the liquid from the sides of the tube and place on the magnet right away until the solution is clear (approx. 1–2 min).
- c. Collect the fragmented mRNA by transferring 10 µL of the supernatant to a nuclease-free 0.2 mL PCR tube.
- d. Add 1 µL QIAseq FastSelect –Globin. Mix thoroughly by pipetting up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes.
- e. Incubate in a thermal cycler with a heated lid as described in Table 14.

Table 14. FastSelect hybridization protocol

Step	Time and temperature
1	2 min at 75°C
2	2 min at 70°C
3	2 min at 65°C
4	2 min at 60°C
5	2 min at 55°C
6	2 min at 37°C
7	2 min at 25°C
8	Hold at 4°C

4. Refer to the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* and immediately proceed to “First Strand cDNA Synthesis.”

Note: “First Strand cDNA Synthesis” is chapter 1.3 in Version 3.1 of the instruction manual.

5. Follow the *NEBNext Ultra II Directional RNA Library Prep Kit for Illumina Instruction Manual* to perform all remaining library construction steps.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –rRNA and/or –Globin with the KAPA RNA HyperPrep Kit

Important points before starting

- The KAPA RNA HyperPrep Kit (Roche cat. no. KK8540 and KK8541) is required for use with this protocol.
- This protocol has been tested with 25 ng to 1 µg of total RNA.
- **Important:** It is highly recommended to dilute the KAPA adapters 1.5-fold compared to what is suggested in the default KAPA protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *KAPA RNA HyperPrep Kit Technical Data Sheet* (KR1350 – v2.17).

Procedure

1. Thaw the tube(s) from the appropriate QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.

2. From the KAPA RNA HyperPrep Kit, prepare the fragmentation and priming mix described in Table 15 at room temperature in a nuclease-free tube.

Table 15. KAPA RNA HyperPrep fragmentation and priming mix

Component	Volume/reaction (μL)
Total RNA (25 ng–1 μg)	9*
Fragment, prime, and elute buffer (2X) [†]	10
Total volume	19

* Reduce volume to 8 μL if removing rRNA and globin.

[†] From KAPA RNA HyperPrep Kit.

3. To the assembled fragmentation and priming mix, add QIAseq FastSelect as follows:
 - **Option 1 (remove rRNA):** Add 1 μ L of the appropriate QIAseq FastSelect –rRNA
 - **Option 2 (remove globin):** Add 1 μ L of QIAseq FastSelect –Globin
 - **Option 3 (remove rRNA and globin):** Add 1 μ L of QIAseq FastSelect –rRNA HMR and 1 μ L of QIAseq FastSelect –Globin
4. Mix thoroughly by gently pipetting the reaction up and down several times and then briefly centrifuge to collect residual liquid from the sides of the tubes

- Incubate in a thermal cycler with a heated lid as described in Table 16 according to your input RNA quality.

Important: Regardless of time and temperature chosen in step 1, steps 2–9 must be performed.

Table 16. Combined KAPA RNA HyperPrep fragmentation and FastSelect hybridization protocol

Input RNA type	Step	Time and temperature
Intact	1 *	Choose: 8 min at 94°C or 6 min 94°C or 6 min at 85°C
Partially degraded	1 †	1–6 min at 85°C
Degraded (e.g., FFPE)	1 ‡	No fragmentation
Steps 2–9 are performed regardless of input RNA quality. They need to be performed whether the RNA is high quality, moderate quality, FFPE, or degraded.	2	2 min at 75°C
	3	2 min at 70°C
	4	2 min at 65°C
	5	2 min at 60°C
	6	2 min at 55°C
	7	2 min at 37°C
	8	2 min at 25°C
	9	Hold at 4°C

* Choose one option, depending if you want a desired mean library insert size of 100–200 bp (8 min at 94°C), 200–300 bp (6 min 94°C), or 300–400 bp (6 min at 85°C).
† For a desired mean library insert size of 100–300 bp.
‡ For a desired mean library insert size of 100–200 bp.

- Refer to the *KAPA RNA HyperPrep Kit Technical Data Sheet* and immediately proceed to “1st Strand Synthesis”, section 3 in v2.17.

7. Follow the *KAPA RNA HyperPrep Kit Technical Data Sheet* to perform all remaining library construction steps.

IMPORTANT



It is highly recommended to dilute the KAPA adapters 1.5-fold compared to what is suggested in the default KAPA protocol.

IMPORTANT



When removing globin, two additional cycles of library amplification need to be performed.

Protocol: FastSelect –Globin with the KAPA mRNA HyperPrep Kit

Important points before starting

- The KAPA mRNA HyperPrep Kit (Roche cat. no. KK8580 and KK8581) is required for use with this protocol.
- **Important:** It is highly recommended to dilute the KAPA adapters 1.5-fold compared to what is suggested in the default KAPA protocol.
- **Important:** When removing globin, two additional cycles of library amplification need to be performed.
- Refer to the *KAPA mRNA HyperPrep Kit Technical Data Sheet* (KR1352 – v5.17).

Procedure

1. Follow the *KAPA mRNA HyperPrep Kit Technical Data Sheet*, “Library Construction Protocol”, section 1 (Reagent Preparation).
2. Follow the *KAPA mRNA HyperPrep Kit Technical Data Sheet*, “Library Construction Protocol”, section 2 (mRNA Capture).
3. Follow the *KAPA mRNA HyperPrep Kit Technical Data Sheet*, “Library Construction Protocol”, section 3 (mRNA Elution, Fragmentation and Priming), steps 3.1–3.4.
4. Thaw the FastSelect –Globin tube from the QIAseq FastSelect kit. Mix by vortexing and then briefly centrifuge to collect residual liquid from the sides of the tubes.
5. In place of step 3.5 in the *KAPA mRNA HyperPrep Kit Technical Data Sheet*, carefully transfer 19 µL of the supernatant containing the eluted, fragmented, and primed RNA into a separate plate or tube.

- To the supernatant, add 1 μ L of QIAseq FastSelect –Globin.
- Mix thoroughly by gently pipetting the reaction up and down several times, and then briefly centrifuge to collect residual liquid from the sides of the tubes.
- Incubate in a thermal cycler with a heated lid as described in Table 17, according to your input RNA quality.

Table 17. FastSelect hybridization protocol

Input RNA type	Step	Time and temperature
Hybridization	2	2 min at 75°C
	3	2 min at 70°C
	4	2 min at 65°C
	5	2 min at 60°C
	6	2 min at 55°C
	7	2 min at 37°C
	8	2 min at 25°C
	9	Hold at 4°C

- Refer to the *KAPA mRNA HyperPrep Kit Technical Data Sheet* and immediately perform step 3.6 (“Place the plate/tube[s] on ice and proceed immediately to 1st Strand Synthesis [step 4]”).

Note: Step 3.6 is found in section 3 in KR1352 – v5.17.

- Follow the *KAPA mRNA HyperPrep Kit Technical Data Sheet* to perform all remaining library construction steps.

IMPORTANT

It is highly recommended to dilute the KAPA adapters 1.5-fold compared to what is suggested in the default KAPA protocol.

IMPORTANT

When removing globin, two additional cycles of library amplification need to be performed.

Troubleshooting Guide

This troubleshooting guide may be helpful in solving any problems that may arise. For more information, see also the Frequently Asked Questions page in our Technical Support Center: www.qiagen.com/FAQ/FAQList.aspx. The scientists in QIAGEN Technical Services are always happy to answer any questions you may have about either the information or protocols in this handbook (for contact information, visit support.qiagen.com).

Comments and suggestions

Adapter-dimer observed in final library QC

Depending on the library kit and RNA input amount, adapter-dimers may be observed.

Perform a second cleanup of the final library using the same beads-to-sample ratio as in the first cleanup.

Appendix A: Species Compatibility

QIAseq FastSelect –rRNA HMR and –Globin species compatibility

In addition to human, mouse, and rat samples, QIAseq FastSelect –rRNA HMR has been tested with a variety of other RNA samples (Table 18). QIAseq FastSelect –Globin has been tested with whole-blood total RNA samples from human, mouse, and rat, wherein 95–99% removal of globin mRNA is routinely observed. Similar to FastSelect –rRNA HMR, FastSelect –Globin may work on species beyond human, mouse, and rat. Importantly, when working with total RNA blood samples, QIAseq FastSelect –rRNA HMR must be used in combination with FastSelect –Globin. When working with mRNA-enriched samples, only FastSelect –Globin needs to be used.

Table 18. Summary of rRNA removal using QIAseq FastSelect –rRNA HMR

Species*	% rRNA reads (no FastSelect)	% rRNA reads (with FastSelect)	% knockdown
Human	91.8	1.6	98
Mouse	92.9	2.2	98
Rat	93.0	2.6	97
Chicken	88.3	16.3	82
Cow	70.0	1.2	98
Cynomolgus monkey (<i>Macaca fascicularis</i>)	90.4	10.6	88
Dog	87.8	9.1	90
Hamster	77.5	3.5	95
Horse	74.1	1.8	98
Pig	87.6	17.3	80
Rabbit	88.8	16.5	81
Sheep	87.3	4.0	95

* We do not recommend using QIAseq FastSelect –rRNA HMR for rRNA removal from roundworm (*Caenorhabditis elegans*) or zebrafish (*Danio rerio*). Laboratory tests suggest minimal rRNA removal from these species using FastSelect –rRNA HMR.

QIAseq FastSelect –rRNA Plant: Types of rRNAs that are removed and species

QIAseq FastSelect –rRNA Plant has been designed to remove cytoplasmic (5.8S, 18S, and 25S), mitochondrial (5S, 18S, and 26S), and chloroplast (4.5S, 5S, 16S, and 23S) rRNA from *Arabidopsis thaliana* and *Arabidopsis lyrata*. In addition to *Arabidopsis* total RNA, QIAseq FastSelect –rRNA Plant has been tested with a variety of other RNA samples (Table 19). Depending on rRNA sequence homology, FastSelect –rRNA Plant will work on additional plant species.

Table 19. Summary of rRNA removal using QIAseq FastSelect –rRNA Plant

Species	% rRNA reads without FastSelect	% rRNA reads with FastSelect	% knockdown
Arabidopsis	96.2	0.8	99
Barley	96.0	7.9	92
Corn	94.7	9.8	90
Cotton	94.1	1.3	99
Flaxseed	95.9	12.6	87
Maple	93.5	10.5	89
Oat	94.8	0.7	99
Potato	95.1	6.5	93
Rice	94.7	8.1	91
Rye	96.2	0.8	99
Sorghum	96.2	20.2	79
Soybean	96.0	6.3	93
Wheat	95.8	5.5	94

QIAseq FastSelect –Mitochondrial Genes Kit

QIAseq FastSelect –Mitochondrial Genes has been designed to remove human, mouse, and rat mitochondrial genes — Human (hg-38): MT-RNR1, MT-RNR2, MT-ND1, MT-ND2, MT-CO1, MT-CO2, MT-ATP8, MT-ATP6, MT-CO3, MT-ND3, MT-ND4L, MT-ND4, MT-ND5, MT-ND6, MT-CYB; Mouse (mm10): mt-Rnr1, mt-Rnr2, mt-Nd1, mt-Nd2, mt-Co1, mt-Co2, mt-Atp8, mt-Atp6, mt-Co3, mt-Nd3, mt-Nd4l, mt-Nd4, mt-Nd5, mt-Nd6, mt-Cytb; Rat (rn6): AY172581.9, AY172581.24, Mt-nd1, Mt-nd2, Mt-co1, Mt-co2, Mt-atp8, Mt-atp6, Mt-co3, Mt-nd3, Mt-nd4l, Mt-nd4, Mt-nd5, Mt-nd6, Mt-cyb.

Species	% mitochondrial genes reads without FastSelect –Mitochondrial Genes Kit	% mitochondrial genes reads with FastSelect –Mitochondrial Genes Kit	% knockdown
Human	5.2	0.1	99
Mouse	7.5	0.5	99
Rat	7.0	0.1	99

Appendix B: FastSelect –rRNA, –Globin mRNA, and Mitochondrial Genes with the QIAseq Stranded RNA Lib Kit

Important points before starting

- This protocol is for combining three different FastSelect panels together in one sample and includes an extra bead cleanup step after RNA fragmentation and before cDNA synthesis.
- The QIAseq Stranded Total RNA Lib Kit (cat. no. 180743 or 180745) is required for use with this protocol.
- This protocol has been tested with 100 ng to 1 µg of total RNA.
- When removing globin, two additional cycles of Hi-Fi library amplification need to be performed.
- Refer to the *QIAseq Stranded RNA Library Kit Handbook* available at www.qiagen.com/HB-2465

Procedure

1. Thaw total RNA on ice. Gently mix, briefly centrifuge to collect residual liquid from the sides of the tubes, and return to ice.
2. Prepare the reagents required for the RNA fragmentation and the appropriate QIAseq FastSelect rRNA, globin, and mitochondrial gene removal.
 - a. Bring FastSelect HMR, FastSelect –Globin, and FastSelect –Mitochondrial Genes Kits and Nuclease-free Water to room temperature.
 - b. Mix by vortexing and then briefly centrifuge.

3. On ice, prepare the fragmentation/depletion reaction according to Table 20. Briefly centrifuge, mix by pipetting up and down 10 times, and centrifuge briefly again.

Note: If setting up more than one reaction, prepare a volume of master mix that is 10% greater than what is required for the total number of reactions.

Table 20. Reaction setup for rRNA + Globin + Mitochondrial Genes depletion

Component	Volume/reaction
Total RNA (100 ng to 1 µg)	Variable
RT Buffer, 5×*	1.5 µL
QIAseq FastSelect HMR	1 µL
QIAseq FastSelect –Globin	1 µL
QIAseq FastSelect –Mitochondrial Genes	1 µL
Nuclease-free water	Bring total reaction volume to 15 µL
Total volume	15 µL*

* From QIAseq Stranded Total RNA Lib Kit.

4. Incubate as described in Table 21, according to input RNA quality and desired insert size.

Important: Regardless of the time and temperature chosen in step 1, steps 2–9 must be performed.

Table 21. Combined QIAseq fragmentation and FastSelect hybridization protocol

Input RNA quality	Step	Insert size approx. 150–250 bp	Insert size approx. 350 bp
High quality (RIN >9)	1 *	15 min at 95°C	3 min at 95°C
Moderate quality (RIN 5–6)	1 *	10 min at 95°C	3 min at 95°C
FFPE or degraded sample (RIN <3)	1 *	No fragmentation†	No fragmentation†
Steps 2–9 are performed regardless of input RNA quality. They need to be performed whether the RNA is high quality, moderate quality, FFPE, or degraded.	2	2 min at 75°C	2 min at 75°C
	3	2 min at 70°C	2 min at 70°C
	4	2 min at 65°C	2 min at 65°C
	5	2 min at 60°C	2 min at 60°C
	6	2 min at 55°C	2 min at 55°C
	7	2 min at 37°C	2 min at 37°C
	8	2 min at 25°C	2 min at 25°C
	9	Hold at 4°C	Hold at 4°C

* Choose one option for the time on step 1 according to the input RNA quality and desired insert size.

† Also suitable for exosomal RNA or RNA of other origin with a size between 80 and 500 bp.

5. Add 19.5 µL (1.3× volume) QIAseq Beads to the 15 µL reaction. Mix thoroughly by vortexing, and then incubate for 5 min at room temperature.
6. Centrifuge in a table-top centrifuge until the beads are completely pelleted (approx. 2 min).

Note: QIAseq Beads should be brought to room temperature and thoroughly mixed by vortexing until a homogenous solution is achieved.

7. Place the tubes/plate on a magnetic rack for 2 min. Once the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.

Note: You can completely avoid transferring beads by pipetting very slowly.

Important: Do not discard the beads, because they contain the RNA of interest.

8. Add 15 μL of Nuclease-free Water and 19.5 μL of QIAseq Bead Binding Buffer (prewarmed to room temperature and mixed by vortexing). Mix vigorously by vortexing, and then incubate for 5 min at room temperature.
9. Centrifuge in a table-top centrifuge until the beads are completely pelleted (approx. 2 min).
10. Place the tubes/plate on a magnetic rack for 2 min. Once the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.

Note: You can completely avoid transferring beads by pipetting very slowly.

Important: Do not discard the beads, because they contain the RNA of interest.

11. With the beads still on the magnetic stand, add 200 μL of 80% ethanol. Wait for 30 s. Carefully remove and discard the wash.
12. Repeat the ethanol wash.

Important: Completely remove all traces of ethanol after this second wash. Remove the ethanol with a 200 μL pipette tip first, and then use a 10 μL pipette tip to remove any residual ethanol that will settle.

13. With the beads still on the magnetic stand, air-dry at room temperature for 6–10 min until all liquid has evaporated but without overdrying the beads.

Note: Visually inspect the pellet to confirm that it is completely dry.

14. Remove the beads from the magnetic stand, and elute the RNA from the beads by adding 31 μL Nuclease-free Water. Mix well by pipetting or vortexing, and then allow the tubes to sit at room temperature for 2 min to appropriately hydrate the beads.
15. Return the tubes/plate to the magnetic rack until the solution has cleared.

- 16. Transfer 29 μ L of the supernatant to clean tubes/plate. Alternatively, the samples can be stored at -90°C to -65°C in a constant-temperature freezer.
- 17. Set up the first-strand synthesis associated with the QIAseq Stranded RNA Lib Kit as described in Table 22. Briefly centrifuge, mix by pipetting up and down 10 times, and then centrifuge briefly again.

Table 22. QIAseq Total RNA Lib Kit first-strand synthesis setup

Component	Volume/reaction (μ L)
RNA from bead cleanup reaction	29
RT Buffer, 5x*	8
Dilute DTT (0.4M)*	1
RT Enzyme*	1
RNase Inhibitor*	1
Total volume	40

* All designated components are from the QIAseq Stranded Total RNA Lib Kit.

- 18. Refer to the *QIAseq Stranded RNA Library Kit Handbook* and immediately proceed to “Protocol: First-strand Synthesis”.
- 19. Follow the *QIAseq Stranded RNA Library Kit Handbook* to perform all remaining library construction steps.

IMPORTANT



When removing globin, HMR rRNA, and mitochondrial genes, two additional cycles of Hi-Fi library amplification need to be performed.

Ordering Information

Product	Contents	Cat. no.
QIAseq FastSelect –rRNA HMR Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports human, mouse, and rat	334386
QIAseq FastSelect –rRNA HMR Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports human, mouse, and rat	334387
QIAseq FastSelect –rRNA HMR Kit (384)	For 384 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports human, mouse, and rat	334388
QIAseq FastSelect –Mitochondrial Genes Kit (24)	For 24 reactions: removal of mitochondrial protein coding genes for human, mouse & rat samples	333282
QIAseq FastSelect –Mitochondrial Genes Kit (96)	For 96 reactions: removal of mitochondrial protein coding genes for human, mouse & rat samples	333286
QIAseq FastSelect Epidemiology Kit	Human/mouse/rat cytoplasmic and mitochondrial rRNA removal reagent, bacteria rRNA removal reagent, fragmentation/hybridization reagent, QIAseq Beads; available in 24, 96, or 384 reactions	333272 333275 333277
QIAseq FastSelect –Globin Kit (24)	For 24 reactions: globin mRNA removal reagent; supports human, mouse, and rat	334376
QIAseq FastSelect –Globin Kit (96)	For 96 reactions: globin mRNA removal reagent; supports human, mouse, and rat	334377
QIAseq FastSelect –Globin Kit (384)	For 384 reactions: globin mRNA removal reagent; supports human, mouse, and rat	334378
QIAseq FastSelect –rRNA/Globin Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent and globin mRNA removal reagent; supports human, mouse, and rat	335376
QIAseq FastSelect –rRNA/Globin Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent and globin mRNA removal reagent; supports human, mouse, and rat	335377
QIAseq FastSelect –rRNA/Globin Kit (384)	For 384 reactions: cytoplasmic and mitochondrial rRNA removal reagent and globin mRNA removal reagent; supports human, mouse, and rat	335378

Product	Contents	Cat. no.
QIAseq FastSelect –rRNA Plant Kit (24)	For 24 reactions: cytoplasmic, mitochondrial and chloroplast rRNA removal reagent; supports plant	334315
QIAseq FastSelect –rRNA Plant Kit (96)	For 96 reactions: cytoplasmic, mitochondrial and chloroplast rRNA removal reagent; supports plant	334317
QIAseq FastSelect –rRNA Plant Kit (384)	For 384 reactions: cytoplasmic, mitochondrial and chloroplast rRNA removal reagent; supports plant	334319
QIAseq FastSelect –rRNA Yeast Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports yeast	334215
QIAseq FastSelect –rRNA Yeast Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports yeast	334217
QIAseq FastSelect –rRNA Yeast Kit (384)	For 384 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports yeast	334219
QIAseq FastSelect –rRNA Worm Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports worm	333242
QIAseq FastSelect –rRNA Worm Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports worm	333245
QIAseq FastSelect –rRNA Fish Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports fish	333252
QIAseq FastSelect –rRNA Fish Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports fish	333255
QIAseq FastSelect –rRNA Fly Kit (24)	For 24 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports fly	333262
QIAseq FastSelect –rRNA Fly Kit (96)	For 96 reactions: cytoplasmic and mitochondrial rRNA removal reagent; supports fly	333265
QIAseq Low Input RNA Library Kit (24)	Includes RNA library preparation reagents for 24 samples, QIAseq Beads and QIAseq Advanced Analysis for the RNA-seq Analysis Portal.	334202
QIAseq Low Input RNA Library Kit (96)	Includes RNA library preparation reagents for 96 samples, QIAseq Beads and QIAseq Advanced Analysis for the RNA-seq Analysis Portal.	334205

Product	Contents	Cat. no.
QIAseq Stranded RNA Lib Kit UDI (24)	For 24 stranded RNA-seq sequencing library preparation reactions: fragmentation, reverse transcription, second-strand synthesis + end-repair + A-addition, adapter ligation, Hi-Fi library amplification, QIAseq Beads for library cleanups, and unique dual index (UDI) adapters	180450
QIAseq Stranded RNA Lib Kit UDI-A (96)/UDI-B (96)/UDI-C (96)/UDI-D (96)	For 96 stranded RNA-seq sequencing library preparation reactions: fragmentation, reverse transcription, second-strand synthesis + end-repair + A-addition, adapter ligation, Hi-Fi library amplification, QIAseq Beads for library cleanups, and unique dual index (UDI) adapters	180451 180452 180453 180454
QIAseq Stranded mRNA Lib Kit UDI (24)	For 24 stranded RNA-seq sequencing library preparation reactions: mRNA enrichment of total RNA, fragmentation, reverse transcription, second-strand synthesis + end-repair + A-addition, adapter ligation, Hi-Fi library amplification, QIAseq Beads for library cleanups, and unique dual index (UDI) adapters	180440
QIAseq Stranded mRNA Lib Kit UDI-A (96)/UDI-B (96)/UDI-C (96)/UDI-D (96)	For 96 Stranded RNA-seq sequencing library preparation reactions: mRNA enrichment of total RNA, fragmentation, reverse transcription, second-strand synthesis + end-repair + A-addition, adapter ligation, Hi-Fi library amplification, QIAseq Beads for library cleanups, and unique dual index (UDI) adapters	180441 180442 180443 180445

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Service or your local distributor.

Document Revision History

Date	Description
June 2021	Initial release
December 2021	Updated the Ordering Information section: Added the QIAseq FastSelect Epidemiology Kit.
January 2026	Updated branding. Added ordering information on QIAseq FastSelect –Mitochondrial Genes Kit. Updated RNA library kit compatibility. Updated links to handbooks.

Limited License Agreement for QIAseq FastSelect products

Use of this product signifies the agreement of any purchaser or user of the product to the following terms:

1. The product may be used solely in accordance with the protocols provided with the product and this Instructions for Use and for use with components contained in the panel only. QIAGEN grants no license under any of its intellectual property to use or incorporate the enclosed components of this panel with any components not included within this panel except as described in the protocols provided with the product, this Instructions for Use, and additional protocols available at www.qiagen.com. Some of these additional protocols have been provided by QIAGEN users for QIAGEN users. These protocols have not been thoroughly tested or optimized by QIAGEN. QIAGEN neither guarantees them nor warrants that they do not infringe the rights of third-parties.
2. Other than expressly stated licenses, QIAGEN makes no warranty that this panel and/or its use(s) do not infringe the rights of third-parties.
3. This panel and its components are licensed for one-time use and may not be reused, refurbished, or resold.
4. QIAGEN specifically disclaims any other licenses, expressed or implied other than those expressly stated.
5. The purchaser and user of the panel agree not to take or permit anyone else to take any steps that could lead to or facilitate any acts prohibited above. QIAGEN may enforce the prohibitions of this Limited License Agreement in any Court, and shall recover all its investigative and Court costs, including attorney fees, in any action to enforce this Limited License Agreement or any of its intellectual property rights relating to the panel and/or its components.

For updated license terms, see www.qiagen.com.

Trademarks: QIAGEN®, Sample to Insight®, QIAseq®, CleanStart®, FastSelect® (QIAGEN Group); Bioanalyzer®, Fragment Analyzer™, TapeStation® (Agilent Technologies, Inc.); Illumina®, TruSeq® (Illumina, Inc.); NEB®, NEBNext®, Ultra™ (New England Biolabs); KAPA®, Roche® (Roche Group); Takara Bio® (Takara Bio Inc.); Watchmaker® (Watchmaker). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

01/2026 HB-2929-003 © 2026 QIAGEN, all rights reserved.

This page intentionally left blank

