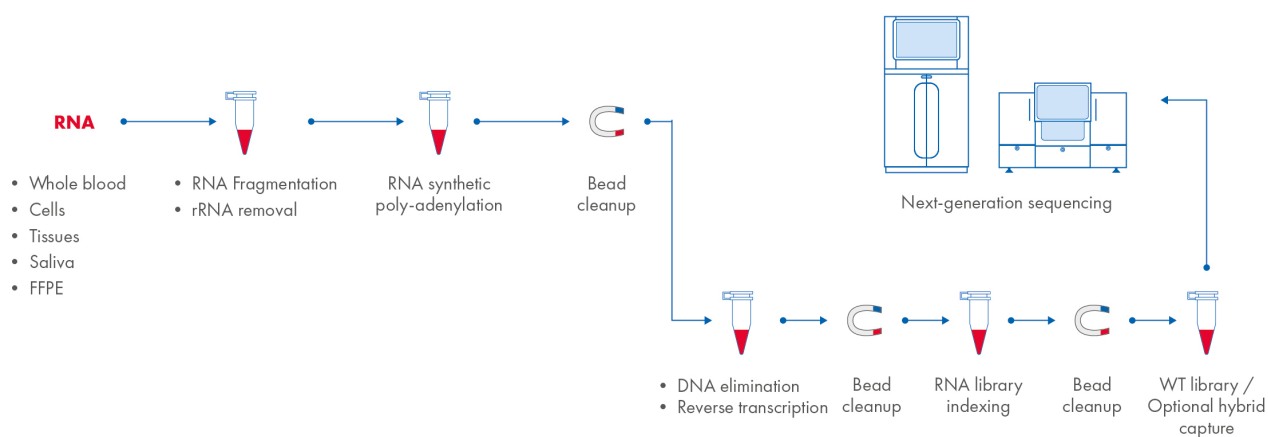


Supplementary Protocol

QIAseq® Multimodal RNA Kit

For the construction of WTS libraries using the QIAseq Multimodal RNA Kit (cat. nos. 335012 and 335015) and QIAseq Multimodal RNA UDI (cat. nos. 334822 and 334825).

QIAseq Multimodal DNA/RNA Library Kit workflow



Equipment and Reagents to be Supplied by User

When working with chemicals, always wear a suitable lab coat, disposable gloves, and protective goggles. For more information, consult the appropriate safety data sheets (SDSs) available from the product supplier.

The following are required:

- 80% ethanol (made fresh daily)*
- Nuclease-free pipette tips and tubes
- 1.5 mL LoBind® tubes (Eppendorf®, cat. no. 022431021)
- PCR tubes (0.2 mL individual tubes [VWR®, cat. no. 20170-012] or tube strips [VWR, cat. no. 93001 118]) or plates
- Ice
- Microcentrifuge
- Thermal cycler
- Magnet for bead cleanups:
 - **Tubes:** MagneSphere® Technology Magnetic Separation Stand (Promega, cat. no. Z5342)
 - **Plates:** DynaMag™-96 Side Magnet (Thermo Fisher Scientific Inc., cat. no. 12331D)
- QIAxcel® Connect System: QIAxcel DNA High Resolution Kit

*Do not use denatured alcohol, which contains other substances, such as methanol or methylethylketone.

Protocol: RNA Fragmentation

Important points before starting

- Set up reactions on ice.
- Do not vortex any reagents or reactions.
- Use MRNA index plates, either MRNA-24X or MRNA-96AX.
- Prepare fresh 80% ethanol daily.
- Ensure that the QIAseq Beads are thoroughly mixed at all times. This necessitates working quickly and resuspending the beads immediately before use. If a delay in the protocol occurs, simply vortex the beads before proceeding.

Procedure

1. Thaw nucleic acid sample(s) on ice. Gently mix, briefly centrifuge to collect residual liquid from the sides of the tubes, and then return to ice.
2. Prepare the reagents required for the fragmentation.
 - a. Thaw 10x FX Buffer and QIAseq FastSelect® –rRNA HMR (8) at room temperature (15–25°C).
 - b. Mix by flicking the tube, and then centrifuge briefly.
3. Dilute an aliquot of QIAseq FastSelect –rRNA HMR 1:10 with Nuclease-free Water. Mix by flicking the tube, and then centrifuge briefly.
4. On ice, prepare the fragmentation mix according to Table 1. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.

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

Table 1. Reaction mix for nucleic acid fragmentation of RNA samples

Component	Volume/reaction (µL)
RNA sample	Variable
10x FX Buffer	2
QIAseq FastSelect –rRNA HMR (diluted 1:10)	1
Nuclease-free Water	17 – Variable
Total	20

5. Program the thermal cycler according to Table 2 and Table 3. Use the instrument's heated lid.

Important: The thermal cycler must be prechilled and paused at 4°C.

Table 2. Incubation conditions for standard RNA fragmentation

Step	Incubation temperature (°C)	Incubation time for RNA (min)
1	4	1
2	95	10
3	75	2
4	70	2
5	65	2
6	60	2
7	55	2
8	37	2
9	25	2
10	4	Hold

Table 3. Incubation conditions for FFPE RNA fragmentation

Step	Incubation temperature (°C)	Incubation time for RNA (min)
1	4	1
2	72	30
3	65	2
4	60	2
5	55	2
6	37	2
7	25	2
8	4	Hold

6. Transfer the tubes/plate wells prepared in step 3 to the prechilled thermal cycler and resume the program.
7. Upon completion, allow the thermal cycler to return to 4°C.
8. Place the samples on ice and immediately proceed to "Protocol: RNA Polyadenylation".

Protocol: RNA Polyadenylation

Important points before starting

- The product from “Protocol: RNA Fragmentation” is the starting material for this protocol.
- Set up reactions on ice.
- Do not vortex any reagents or reactions.

Procedure

1. Prepare the reagents required for the polyadenylation.
 - a. Thaw PAP Dilution Buffer, 10x, and ATP Solution on ice.
 - b. Mix by flicking the tube, and then centrifuge briefly.

Note: Just before use, remove T4 Polynucleotide Kinase and PAP Enzyme from the freezer, and place them on ice. After use, immediately return the enzymes to the freezer.
2. Dilute PAP Enzyme from 5 U/μL to 1 or 2 U/μL as follows:
 - a. Prepare 1x PAP Dilution Buffer by diluting 2 μL of the 10x PAP Dilution Buffer with 18 μL Nuclease-Free Water.
 - b. Use the 1x PAP Dilution Buffer to dilute an aliquot of the PAP Enzyme from 5 U/μL to 1 or 2 U/μL. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.
 - c. For RNA input ≥80 ng, use 2 U/μL.
3. Prepare the RNA polyadenylation mix according to Table 4. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.

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

Table 4. Reaction mix for RNA polyadenylation

Component	Volume/reaction (μL)
Fragmentation reaction (already in tube)	20
ATP Solution	1.25
T4 Polynucleotide Kinase	1
Diluted PAP Enzyme (1 or 2 U/μL)*	1
Nuclease-Free Water	1.75
Total	25

*Ensure PAP Enzyme has been diluted from its stock 5 U/μL concentration to 1 or 2 U/μL using 1x PAP Dilution Buffer. For RNA input ≥80 ng, use 2 U/μL.

4. Incubate the reactions in a thermal cycler according to Table 5. Use the instrument's heated lid.

Table 5. Incubation conditions for RNA polyadenylation

Step	Temperature (°C)	Time (min)
1	4	1
2	30	10
3	4	Hold

5. After the reaction, add 55 μL of Nuclease-Free Water to make a total volume of 80 μL .
6. Add 104 μL QIAseq Beads and mix by vortexing.
7. Incubate for 5 min at room temperature.
8. Place the tubes/plate wells on a magnetic rack for 2 min (for tubes) or 10 min (for plates). Once the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.

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

Optional: For plates, the following may improve performance:

- a. After 8 min, remove 100 μL supernatant. Leave it on the magnetic stand for 2 min and remove 100 μL supernatant.
- b. Briefly centrifuge the plate, return it to the magnetic stand for 1 min, and use a 10 μL pipette to remove the remaining supernatant.
- c. Spin down briefly and remove the residual.
9. Resuspend in 40 μL of Nuclease-Free Water and add 52 μL QIAseq Beads. Mix well and incubate for 5 min at room temperature.
10. With the beads still on the magnetic stand, add 200 μL of 80% ethanol. Rotate the tube 2–3 times or move the plate side to side between the 2 column positions of the magnet to wash the beads. Carefully remove and discard the wash.
11. Repeat the ethanol wash.

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

12. With the beads still on the magnetic stand, air-dry at room temperature for 5–10 min.

Note: Visually inspect the pellet to confirm that it is completely dry. If the pellet is not completely dry, additional drying time may be required.

13. Remove the beads from the magnetic stand, and then elute the RNA from the beads by adding 12 μL of Nuclease-Free Water. Mix well by pipetting.
14. Return the tubes/plate wells to the magnetic rack until the solution has cleared.
15. Transfer 8.4 μL of the eluate to clean tubes/plate wells.
16. Proceed to “Protocol: RNA Reverse Transcription”.

Alternatively, the samples can be stored at -30°C to -15°C in a constant-temperature freezer.

Protocol: RNA Reverse Transcription

Important points before starting

- The product from “Protocol: RNA Polyadenylation” is the starting material for this protocol.
- Set up reactions on ice.
- Do not vortex any reagents or reactions.
- Prepare fresh 80% ethanol daily.
- Ensure that the QIAseq Beads are thoroughly mixed at all times. This necessitates working quickly and resuspending the beads immediately before use. If a delay in the protocol occurs, simply vortex the beads before proceeding.

Procedure

1. Prepare the reagents required for reverse transcription.
 - a. Thaw US RT Buffer, 5×; DTT (100 mM); dNTP (10 mM); MM RNA RT Primer; Multimodal RT Enhancer; and MM RNA TSO at room temperature.
 - b. Mix by flicking the tubes and then centrifuge briefly.

Note: Just before use, remove the RNase Inhibitor and EZ Reverse Transcriptase from the freezer and place them on ice. After use, immediately return the enzymes to the freezer.

2. To the treated sample, prepare the reverse transcription mix according to Table 6. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.

Note: If setting up more than 1 reaction, prepare a volume of master mix 10% greater than what is required for the total number of reactions.

Table 6. Reaction mix for reverse transcription

Component	Volume/reaction (μL)
Treated Sample (from previous section)	8.4
US RT Buffer, 5×	4
DTT (100 mM)	0.5
dNTP (10 mM)	2
MM RNA RT Primer	1.6
MM RNA TSO	1.6
Multimodal RT Enhancer	0.4
RNase Inhibitor	0.5
EZ Reverse Transcriptase	1
Total	20

3. Incubate the reactions in a thermal cycler according to Table 7. Use the instrument’s heated lid.

Table 7. Incubation conditions for reverse transcription

Step	Incubation temperature (°C)	Incubation time (min)
1	4	1
2	25	5
3	42	90
4	70	10
5	4	1
6	4	Hold

4. Add 20 µL Nuclease-Free Water and 44 µL QIAseq Beads and mix by vortexing or by pipetting up and down several times.
5. Incubate for 5 min at room temperature.
6. Place the tubes/plate wells on a magnetic rack for 2 min (for tubes) or 10 min (for plates).
After the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.
Important: Do not discard the beads because they contain the DNA of interest.
7. With the beads still on the magnetic stand, add 200 µL of 80% ethanol. Rotate the tube 2–3 times or move the plate side to side between the 2 column positions of the magnet to wash the beads. Carefully remove and discard the wash.
8. Repeat the ethanol wash.
Important: Completely remove all traces of the ethanol after this second wash. Remove the ethanol with a 200 µL pipette tip first, spin down briefly, and then use a 10 µL pipette tip to remove any residual ethanol.
9. With the beads still on the magnetic stand, air-dry at room temperature for 5–10 min.
Note: Visually inspect the pellet to confirm that it is completely dry. If the pellet is not completely dry, additional drying time may be required. Ethanol carryover to the next universal PCR step will affect PCR efficiency.
10. Remove the beads from the magnetic stand, and then elute the DNA from the beads by adding 42 µL Nuclease-Free Water.
11. Return the tubes/plate wells to the magnetic rack until the solution has cleared.
12. Transfer 40 µL of the eluate to clean tubes/plate wells.
13. Add 44 µL QIAseq Beads and mix by vortexing or by pipetting up and down several times.
14. Incubate for 5 min at room temperature.
15. Place the tubes/plate wells on a magnetic rack for 2 min (for tubes) or 10 min (for plates). After the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.
Important: Do not discard the beads because they contain the DNA of interest.
16. With the beads still on the magnetic stand, add 200 µL of 80% ethanol. Rotate the tube 2–3 times or move the plate side to side between the 2 column positions of the magnet to wash the beads. Carefully remove and discard the wash.
17. Repeat the ethanol wash.

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

18. With the beads still on the magnetic stand, air-dry at room temperature for 5–10 min.

Note: Visually inspect the pellet to confirm that it is completely dry. If the pellet is not completely dry, additional drying time may be required. Ethanol carryover to the next universal PCR step will affect PCR efficiency.

19. Remove the beads from the magnetic stand, and then elute the DNA from the beads by adding 23 µL Nuclease-Free Water.
20. Return the tubes/plate wells to the magnetic rack until the solution has cleared.
21. Transfer 20 µL of the eluate to clean tubes/plate wells.
22. Proceed to “Protocol: RNA Library Indexing”.

Alternatively, the samples can be stored at –30°C to –15°C in a constant-temperature freezer.

Protocol: RNA Library Indexing

Important points before starting

- The product from "Protocol: RNA Reverse Transcription" is the starting material for this protocol.
- Set up reactions on ice.
- Do not vortex any reagents or reactions.
- Use MRNA index plates, either MRNA-24X or MRNA-96AX.
- Prepare fresh 80% ethanol daily.
- Ensure that the QIAseq Beads are thoroughly mixed at all times. This necessitates working quickly and resuspending the beads immediately before use. If a delay in the protocol occurs, simply vortex the beads before proceeding.

Procedure

1. Prepare the reagents required for RNA Library indexing.
 - a. Thaw HiFi Ultra Buffer, 5x; MM F-R Primer Mix, and MRNA-24X or MRNA-96AX index plate at room temperature.
 - b. Mix by either flicking the tube or vortexing the index plate, and then centrifuge briefly. Plate should be centrifuged at $1000 \times g$ for 1 min.

Note: Just before use, remove HiFi Ultra Polymerase from the freezer and place on ice. After use, immediately return the enzyme to the freezer.

2. Prepare the reactions according to Table 8. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.

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

Table 8. Reaction mix for RNA indexing

Component	Volume/reaction (µL)
Sample aliquot (from "Protocol: RNA Reverse Transcription")	20
HiFi Ultra Buffer, 5x	8
MM F-R Primer Mix†	1.6
Nuclease-Free Water	3.9
Well from MRNA-24X or MRNA-96AX index plate*	2.5
HiFi Ultra Polymerase	4
Total	40

*Ensure proper technique to prevent cross-contamination. Additionally ensure that every sample has a unique index and that no well is used twice.

† If the input is ≤ 10 ng, reduce the volumes of both the MM F-R Primer Mix and the corresponding well from the MRNA-24X or MRNA-96X index plate to half of the standard amount.

3. Program a thermal cycler as described in Table 9, using cycle numbers described in Table 10.

Important: Set up the ramp rate of the thermal cycler at $\leq 2^{\circ}\text{C}/\text{s}$.

Table 9. Cycling conditions for RNA indexing

Step	Time	Temperature ($^{\circ}\text{C}$)
Hold	2 min	98
2-step cycling		
Denaturation	20 s	98
Annealing	1 min	60
Extension	1 min	72
Cycle number	See Table 10	
Hold	3 min	72
Hold	∞	4

Table 10. Cycle number recommendations for RNA indexing, based on original sample input

RNA Input (ng)	Standard Sample (cycle)	FFPE sample (cycle)
5	25	–
10	23	–
20	22	–
80	20	22
250	19	20
500	–	19

4. After the reaction is complete, add 36 μL QIAseq Beads, and then mix by vortexing or pipetting up and down several times.
5. Incubate for 5 min at room temperature.
6. Place the tubes/plate wells on a magnetic rack for 2 min (for tubes) or 10 min (for plates) to separate the beads from the supernatant.

After the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant.

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

7. With the beads still on the magnetic stand, add 200 μL of 80% ethanol. Rotate the tube 2–3 times or move the plate side to side between the 2 column positions of the magnet to wash the beads. Carefully remove and discard the wash.
8. Repeat the ethanol wash twice, for a total of 3 washes.

Important: Completely remove all traces of the ethanol after the third wash. Remove the ethanol with a 200 μL pipette tip first, spin down briefly, and then use a 10 μL pipette tip to remove any residual ethanol.

9. With the beads still on the magnetic stand, air-dry at room temperature for 5–10 min.

Note: Visually inspect the pellet to confirm that it is completely dry. If the pellet is not completely dry, additional drying time may be required.

10. Remove the beads from the magnetic stand, and then elute the DNA from the beads by adding 24 μ L Nuclease-Free Water. Mix well by pipetting.
11. Return the tube/plate wells to the magnetic rack until the solution has cleared.
12. Transfer 22 μ L of the supernatant to clean tubes/plate wells.
13. The library is now ready for sequencing or hybrid capture. Proceed to "Recommendations: Library QC & Quantification" in *QIAseq Multimodal DNA/RNA Library Kit Handbook* (www.qiagen.com/HB-3579). Alternatively, the samples can be stored at -30°C to -15°C in a constant-temperature freezer.

Document Revision History

Date	Description
06/2025	Initial release
03/2026	In "Protocol: RNA Polyadenylation": modified protocol steps. In "Protocol: RNA Reverse Transcription": modified protocol steps and change in tables for reverse transcription reaction mix and incubation conditions. In "Protocol: RNA Library Indexing": change in tables for programming a thermal cycler.

Limited License Agreement for QIAseq® Multimodal RNA Kit

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®, FastSelect®, QIAseq®, QIAxcel® (QIAGEN Group); VWR® (Avantor®, Inc.); LoBind® (Eppendorf AG); MagneSphere® (Promega Corporation); DynaMag™ (Thermo Fisher Scientific Inc.). Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

03/2026 HB-3705-002 © 2026 QIAGEN, all rights reserved.

This page intentionally left blank.

This page intentionally left blank.

