



June 2026

QIAseq[®] xHYB CGP DNA/RNA Panel Handbook

For library preparation and hybridization capture–based enrichment
for comprehensive genomic profiling from FFPE samples

Table of Contents

Kit Contents	4
Shipping and Storage	9
Intended Use	9
Safety Information	10
Quality Control	10
Introduction	11
Principle and procedure	12
QIAseq One-4-All Blocking Oligos	13
QIAseq N Blocking Oligos	13
QIAseq xHYB CGP Panel	13
Description of protocols	14
Equipment and Reagents to be Supplied by User	15
Important Notes	16
Protocol: Nucleic Acid Fragmentation	18
Protocol: RNA Polyadenylation	21
Protocol: DNA Ligation	24
Protocol: DNA Library Indexing	28
Protocol: RNA Reverse Transcription	32
Protocol: RNA Library Indexing	37
Protocol: Target Enrichment from Indexed Whole-genome Libraries Using QIAseq xHYB	
Human Hybrid Capture Panels	41
Pool libraries and dry down the indexed library pool	41
Hybridization capture	44
Bind hybridized targets to Streptavidin beads	46
Postcapture amplification	50
Recommendations: Library QC and Quantification	54
Sequencing Guidelines	55
Pooling and multiplexing guidelines	55
Sequencing	57

Sequencing setup on Illumina instruments 58

Troubleshooting Guide 59

Ordering Information 61

Document Revision History 63

Kit Contents

QIAseq xHYB CGP DNA Panel	(24)	(96)
Catalog no.	333122	333125
No. of hybridization reactions	3	12
Typical no. of samples	24	96
xHYB Probe panel for DNA target enrichment for comprehensive genomic profiling	12 µL	48 µL
QIAseq xHYB CGP RNA Panel	(24)	(96)
Catalog no.	334122	334125
No. of hybridization reactions	3	12
Typical no. of samples	24	96
xHYB Probe panel for RNA target enrichment for comprehensive genomic profiling	12 µL	48 µL
QIAseq xHYB CGP DNA/RNA Panels	(24)	(96)
Catalog no.	335122	335125
No. of hybridization reactions	3	12
Typical no. of samples	24	96
xHYB Probe panel for DNA and RNA target enrichment for comprehensive genomic profiling	12 µL	48 µL

QIAseq xHYB Human Reagent Kit	(24)	(96)
Catalog no.	333430	333195
No. of hybridization reactions	3	12
Typical no. of samples	24	96

Box 1 of 2		
One-4-All Blocking Oligos	24 µL	96 µL
One-4-All Blocking Solution	15 µL	60 µL
Fast Hybridization Solution	60 µL	240 µL
Vapor-Lock	500 µL	500 µL
Post Hybrid-Capture PCR Mix (2×)	660 µL	660 µL
Primer Mix Illumina Library Amplification 12 rxn	20 µL	3 × 20 µL
RNase-Free Water	1.5 mL	1.5 mL
Box 2 of 2		
Streptavidin Binding Beads	300 µL	1.2 mL
Post Capture Binding Buffer	2 × 1.2 mL	9.6 mL
Wash Buffer A	2 × 675 µL	5.4 mL
Wash Buffer B	2 × 1.05 mL	8.4 mL

QIAseq Multimodal DNA/RNA Library Kit	(24)	(96)
Catalog no.	334842	334845
Number of samples	24	96
10x FX Buffer	48 µL	225 µL
5x WGS FX Mix	96 µL	475 µL
QIAseq FastSelect® –rRNA HMR (8)	12 µL	12 µL
Side Reaction Reducer	48 µL	192 µL
FG Solution	170 µL	170 µL
ATP Solution	36 µL	290 µL
PAP Enzyme	96 µL	96 µL
PAP Dilution Buffer	192 µL	192 µL
T4 Polynucleotide Kinase	125 µL	125 µL
DNA Ligase	2 × 120 µL	1100 µL
UPH Ligation Buffer, 2.5x	1152 µL	3 × 1500 µL
Buffer GE2 gDNA elimination buffer	72 µL	250 µL
US RT Buffer, 5x	150 µL	450 µL
100 mM DTT	20 µL	80 µL
10 mM dNTP	55 µL	235 µL
EZ Reverse Transcriptase	36 µL	2 × 96 µL
RNase Inhibitor	12 µL	96 µL
Nuclease-Free Water	10 mL	10 mL
QIAseq Beads	10 mL	55 mL

QIAseq Multimodal DNA/RNA UDI (24)
(DNA and RNA indexing for 24 samples)
Catalog no.

(24)
334852

MDNA24X (QIAseq Multimodal MDNA Index Plate 24)	1
Each plate allows indexing of 24 DNA samples. Each well in the plate is single-use and contains 9 µL of UDI index pairs for DNA libraries.	
MRNA-24X (QIAseq Multimodal MRNA Index Plate 24)	1
Each plate allows indexing of 24 RNA samples. Each well in the plate is single-use and contains 9 µL of UDI index pairs for RNA libraries.	
Non-UMI Adapter	60 µL
UMI Adapter	60 µL
MM RNA RT Primer	38.4 µL
MM RNA TSO Primer	38.4 µL
HiFi Ultra Buffer	384 µL
HiFi Ultra Polymerase	192 µL
MM F-R Primer Mix	77 µL
Multimodal RT Enhancer	12 µL

QIAseq Multimodal DNA/RNA UDI Set A
(DNA and RNA indexing for 96 samples)
Catalog no.

(96)
334855

MDNA-96AX (QIAseq Multimodal MDNA Index 96 Set A)	1
Each plate allows indexing of 96 DNA samples. Each well in the plate is single-use and contains 9 µL of UDI index pairs for DNA libraries.	
MRNA-96AX (QIAseq Multimodal MRNA Index 96 Set A)	1
Each plate allows indexing of 96 RNA samples. Each well in the plate is single-use and contains 9 µL of UDI index pairs for RNA libraries.	
Non-UMI Adapter	275 µL
UMI Adapter	275 µL
MM RNA RT Primer	184 µL
MM RNA TSO Primer	184 µL
HiFi Ultra Buffer	1536 µL
HiFi Ultra Polymerase	850 µL
MM F-R Primer Mix	340 µL
Multimodal RT Enhancer	48 µL

Product	Description	Cat. no.
QIAseq N Blocking Oligos (12)	Oligos designed to bind and block all Illumina Nextera-compatible NGS adapters for 12 reactions of hybridization capture	334871

Shipping and Storage

QIAseq xHYB CGP Panels are shipped on dry ice. Store spike-in panels immediately upon receipt at -30°C to -15°C in a constant-temperature freezer. The QIAseq xHYB Human Reagent Kit is shipped on a cool pack (4°C). Store the automation kit immediately upon receipt at $2-8^{\circ}\text{C}$.

QIAseq Multimodal DNA/RNA Library Kit (except QIAseq Beads and Nuclease-Free Water) is shipped on dry ice and should be stored immediately upon receipt at -30°C to -15°C in a constant-temperature freezer. QIAseq Beads and Nuclease-Free Water are shipped in a separate box on cold packs and should be stored immediately upon receipt at $2-8^{\circ}\text{C}$.

QIAseq Multimodal DNA/RNA Index Kits are shipped on dry ice and should be stored immediately upon receipt at -30°C to -15°C in a constant-temperature freezer.

Under these conditions, the components are stable until the expiry date indicated on the label.

Intended Use

QIAseq xHYB Panels, QIAseq xHYB Human Reagent Kits, QIAseq Multimodal DNA/RNA Library and Index Kits 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 components.

Quality Control

In accordance with QIAGEN ISO-certified Quality Management System, each lot of QIAseq xHYB CGP DNA/RNA Panels is tested against predetermined specifications to ensure consistent product quality.

Introduction

The QIAseq xHYB CGP Panel represents a significant advancement in comprehensive genomic profiling, tailored to enhance cancer research and diagnostics through its multiomic and multimodal capabilities.

This robust panel targets over 720 genes, ensuring full coverage of coding regions and microsatellite instability (MSI) markers, critical for precise tumor profiling and biomarker identification. Its compatibility with cell-free DNA (cfDNA) facilitates the accurate detection of gene fusions, essential for a deeper understanding of cancer biology. The panel's multiomic capability integrates both DNA and RNA analyses, empowering researchers to explore tumor heterogeneity and gene expression dynamics. Customization options further enhance its versatility, allowing users to add specific targets or utilize an optional RNA panel that focuses on over 230 genes.

The multimodal library preparation streamlines both DNA and RNA workflows, optimizing resource use and increasing efficiency, while flexible hybridization times and automation options improve workflow adaptability for diverse research applications.

In summary, the QIAseq xHYB CGP Panel serves as a powerful tool for advancing oncology through comprehensive multiomic genomic analysis, providing valuable insights that can lead to the discovery of new biomarkers and therapeutic targets in cancer, meeting the growing demand for accessible and effective genomic profiling solutions in the evolving landscape of cancer research.

Principle and procedure

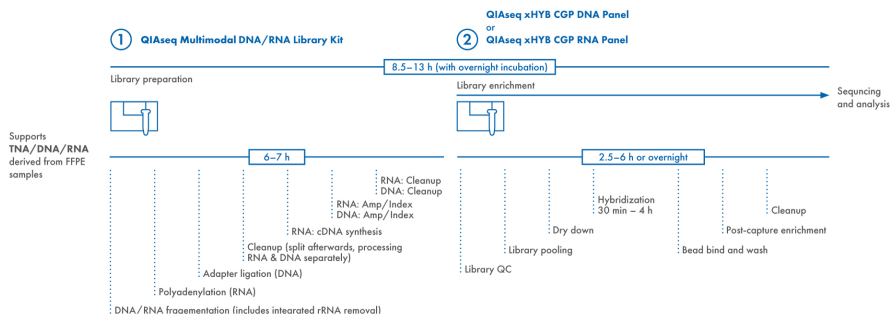


Figure 1. Workflow principle of the QIAseq xHYB CGP Panels. **Step 1:** Generation of indexed libraries using QIAseq Multimodal DNA/RNA Library Kit. **Step 2:** Enrichment of xHYB CGP Targets.

Step 1: Generation of indexed libraries

QIAseq Multimodal DNA/RNA Lib Kit enables construction of WGS and WTS libraries. We recommend using minimum of 40 ng DNA and 80 ng RNA in good quality for CGP DNA/RNA Panels. The following reactions occur in a streamlined, single-tube workflow.

Step 2: Enrichment of human xHYB targets

During xHYB enrichment, indexed library fragments are bound to biotinylated double-stranded DNA capture probes with highly flexible hybridization times ranging from 30 min to overnight incubation. Bound fragments are immobilized on streptavidin beads and non-targeted fragments are washed away. Enriched library fragments are amplified using a proprietary postcapture amplification mix that allows even amplification of DNA regions with vastly different GC contents, which minimizes sequencing bias caused by PCR.

QIAseq One-4-All Blocking Oligos

QIAseq xHYB Human Hybrid Capture Panels include One-4-All Blocking Oligos designed to effectively bind and block all Illumina TruSeq®-compatible NGS adapters with index sequence length of up to 12 nt. All QIAseq Dual-Index Y-Adapters (Unique and Combinatorial) are fully compatible with One-4-All Blocking Oligos for hybridization capture. Using these blocking oligos, the target capture specificity is largely enhanced by preventing capture of library fragments via their adapter sequences during the hybridization reaction.

QIAseq N Blocking Oligos

The RNA libraries generated using the QIAseq Multimodal DNA/RNA Library Kit contain the Nextera® adapter sequence. QIAseq N Blocking Oligos are available (cat. no. 334871, purchased separately) for effectively binding and blocking Nextera-compatible NGS adapters. Using these blocking oligos, the target capture specificity is largely enhanced by preventing capture of library fragments via their adapter sequences during the hybridization reaction.

QIAseq xHYB CGP Panel

QIAseq xHYB CGP DNA panel is a Comprehensive Genomic Profiling assay in with 729 genes enabling thorough assessment of all DNA variant types, including SNVs, CNVs, indels, as well as MSI and TMB. It can be combined with QIAseq xHYB CGP RNA panel targeting 274 genes for fusion detection. This broad and deep analysis supports advanced cancer research applications.

Description of protocols

The QIAseq Multimodal DNA/RNA Library Kit is for use with QIAseq xHYB CGP Panel protocols describes the generation of whole genome and whole transcriptome libraries using an optimized workflow. Alternatively, you may use the QIAseq FX DNA Library Kit, and refer to its respective handbook at www.qiagen.com/QseqFX-DNA

The QIAseq Multimodal DNA/RNA Library Kit can also be used for the generation of DNA-only or RNA-only libraries if needed.

The QIAseq xHYB CGP Panel protocol describes the hybridization capture workflow for any human xHYB panel and is suitable for input libraries generated using QIAseq Multimodal DNA/RNA Library Kit, QIAseq FX DNA Library Kits, QIAseq Ultralow Input Library Kits, or alternative methods.

Equipment and Reagents to be Supplied by User

- QIAseq Beads (cat. nos. 333923, 333903, and 333927) for bead-based library purification
- 100% ethanol (ACS grade)
- Nuclease-Free water
- Buffer EB (cat. no. 19086)
- PCR tubes or plates
- Pipette tips and pipettes
- DNA LoBind tubes (from Axygen™ or Eppendorf®)
- Vortexer
- Microcentrifuge
- Magnetic racks for magnetic beads separation (e.g., Thermo Fisher Scientific DynaMag™ Magnet)
- Thermocycler with heated lid
- Capillary electrophoresis device, e.g., QIAGEN QIAxcel®, Agilent® Bioanalyzer®, or similar to evaluate the DNA fragmentation profile (optional)
- 2 heating blocks for 1.5–2.0 mL tubes
- An evaporation device (e.g., Eppendorf® Concentrator Plus, cat. no. 305000100) equipped with a rotor for 1.5 mL tube and/or a rotor for 0.2 mL PCR tubes/plates.

Important Notes

General precautions

- Use good laboratory practices to minimize cross-contamination of nucleic acid products.
- Always use PCR tubes, microcentrifuge tubes, and pipette tips that are certified sterile and DNase-free and RNase-free.
- Before starting, wipe down the work area and pipettes with an RNase and DNA cleaning product.
- For consistent hybridization capture and amplification, ensure the thermal cycler used in this protocol is in good working condition and has been calibrated within the manufacturer's specifications.
- Read the entire protocol before beginning. Take note of stopping points where samples can be frozen at -30°C to -15°C and plan your workflow accordingly.

Quality control of input libraries

- Control and quantify each indexed library using a capillary electrophoresis method such as using High Sensitivity DNA chip with the QIAxcel Connect System, Agilent 2100 Bioanalyzer, or TapeStation®.
- A typical indexed whole genome library has an average size distribution centered around 400 bp and is free of adapter dimers at a typical yield of >500 ng (refer to Figure 3 in "Recommendations: Library QC and Quantification").
- Libraries from physically sheared DNA will reflect the size distribution of the sheared input DNA.

Indexing recommendations

- Use QIAseq Multimodal DNA/RNA UDI to generate whole genome and/or whole transcriptome libraries for hybridization capture.
- Use different sample indexes for all samples within a hybridization capture pool.
- Consider using different sample indexes also across hybridization capture pools to allow post capture multiplexing depending on your sequencing instrument.

Protocol: Nucleic Acid Fragmentation

Important points before starting

- This protocol describes fragmentation of nucleic acids samples.
- Ensure the nucleic acid sample is in buffer with <0.1 mM EDTA.

Recommended starting amounts of nucleic acid:

- For DNA, use a minimum of 40 ng in good quality with the QIAseq xHYB Kit CGP panel and reagent kits.
 - For RNA, use a minimum of 80 ng in good quality with the QIAseq xHYB Kit CGP panel and reagent kits.
- Set up reactions on ice.
 - Unless specifically indicated, do not vortex any reagents or reactions.

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 10× FX Buffer, QIAseq FastSelect –rRNA HMR (8) at room temperature (15–25°C).
 - b. Mix by flicking the tube, and then centrifuge briefly.

Note: Just before use, remove 5× WGS FX Mix and Side Reaction Reducer from the freezer and place them on ice. After use, immediately return the enzymes to the freezer.

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 FFPE samples

Component	Volume/reaction
DNA (see input recommendation in the “Important points before starting” section)	Variable A
RNA (see input recommendation in the “Important points before starting” section)*	Variable B
10x FX Buffer	2 µL
5x WGS FX Mix	4 µL
QIAseq FastSelect –rRNA HMR (diluted 1:10)	1 µL
Side Reaction Reducer	1.6 µL
Nuclease-Free Water	11.4 µL – Variable A – Variable B
Total	20 µL

*Instead of adding DNA and RNA separately, add total nucleic acid containing both DNA and RNA. If adding total nucleic acid, base the input amount on DNA.

5. Program the thermal cycler according to Table 2. Use the instrument’s heated lid.

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

Table 2. Incubation conditions for nucleic acid fragmentation

Step	Incubation temperature (°C)	Incubation time for FFPE DNA (min)
1	4	1
2	32	12
3	65	30
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: Nucleic Acid Fragmentation and End Repair ” is the starting material for this protocol if input is both DNA and RNA, or RNA-only.
- For DNA-only input, skip this protocol and go directly to “Protocol: DNA Ligation”.
- 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 each reagent by flicking the tube approx. 7 to 8 times, 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 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 2 U/μL. Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.

Dilute PAP Enzyme from 5 U/μL to 1 U/μL as follows:

- a. Prepare 1 × PAP Dilution Buffer by diluting 2 µL of the 10× PAP Dilution Buffer with 18 µL Nuclease-free Water.
 - b. Use the 1 × PAP Dilution Buffer to dilute an aliquot of the PAP Enzyme from 5 U/µL to 1 U/µL (if RNA input <100 ng) or 2 U/µL (if RNA input ≥100 ng). Briefly centrifuge, mix by pipetting up and down 20 times, and then briefly centrifuge again.
3. Prepare the RNA polyadenylation mix according to Table 3Table 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 3. 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 (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 2 U/µL using 1 × PAP Dilution Buffer.

Table 4. Reaction mix for RNA polyadenylation

Component	Volume/reaction (µL)	
	RNA input <100 ng	RNA input ≥100 ng
Fragmentation reaction (already in tube)	20	20

Table 4. Reaction mix for RNA polyadenylation (continued)

Component	Volume/reaction (μL)	
	RNA input <100 ng	RNA input ≥100 ng
ATP Solution	1.25	1.25
T4 Polynucleotide Kinase	1	1
Diluted PAP Enzyme (1 U/μL)*	1	N/A
Diluted PAP Enzyme (2 U/μL)*	N/A	1
Nuclease-Free Water	1.75	1.75
Total	25	25

* Ensure PAP Enzyme has been diluted from its stock 5 U/μL concentration to 1 U/μL or 2 U/μL using 1× PAP Dilution Buffer.

- 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

- Upon completion, place the reactions on ice and proceed to "Protocol: DNA Ligation", next section.
- Upon completion, place the reactions on ice and proceed to
 - "Protocol: DNA Ligation" if input is both DNA and RNA; or
 - "Protocol: Clean up Ligation or Polyadenylation with QIAseq bead" if input is RNA-only.

Protocol: DNA Ligation

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 the DNA ligation.
 - a. Thaw UMI Adapter and UPH Ligation Buffer, 2.5×, at room temperature.
Note: The UMI Adapter contains UMI.
 - b. Mix by flicking the tube, and then centrifuge briefly.
Note: Just before use, remove DNA Ligase from the freezer and place on ice. After use, immediately return the enzyme to the freezer.
2. Prepare the DNA ligation 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 one reaction, prepare a volume of master mix 10% greater than what is required for the total number of reactions.

Table 6. Reaction mix for DNA ligation

Component	Volume/reaction (µL)
RNA polyadenylation reaction (already in tube)	25
UPH Ligation Buffer, 2.5x	40
UMI Adapter*	2.5
DNA Ligase	10
Nuclease-Free Water	22.5
Total	100

*For FFPE samples with input <=100 ng, use UMI Adapter with 1:10 dilution.

3. Incubate the reactions in a thermal cycler according to Table 7.

Important: Do not use the heated lid.

Table 7. Incubation conditions for DNA ligation

Step	Temperature (°C)	Time (min)
1	4	1
2	20	15
3	4	Hold

4. Add 130 µL QIAseq Beads, and then mix by vortexing.
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).

Once the solution has cleared, with the beads still on the magnetic stand, carefully remove and discard the supernatant. Spin down briefly, and then use a 10 µL pipette tip to remove any residual supernatant.

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

Note: For plates, the following may improve performance. After 8 min, remove 90 μ L supernatant. Leave it on the magnetic stand for 2 min and remove 90 μ L supernatant. Then, briefly centrifuge the plate, return it to the magnetic stand for 1 min, and use a 10 μ L pipette to remove the remaining supernatant.

7. Resuspend the beads with 40 μ L Nuclease-Free Water. Then add 52 μ L QIAseq Beads and mix well by pipetting.
8. Incubate for 5 min at room temperature.
9. Place the tubes/plate 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 DNA of interest.

10. With the beads still on the magnetic stand, add 200 μ L 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 μ L pipette tip first, spin down briefly, and then use a 10 μ L 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 DNA from the beads by adding 16 μ L Nuclease-Free Water. Mix well by pipetting.
14. Return the tubes/plate wells to the magnetic rack until the solution has cleared.

15. For each sample, aliquot 6.8 μ L of the eluate into two separate clean tubes/plate wells.

Aliquot 1: Proceed to “Protocol: DNA Library Indexing” on the next page.

Aliquot 2: Proceed to “Protocol: RNA Reverse Transcription” on page 32

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

Protocol: DNA Library Indexing

Important points before starting

- The starting material is a 6.8 μ L sample aliquot from “Protocol: DNA Ligation”.
- Set up reactions on ice.
- Do not vortex any reagents or reactions.
- Use MDNA index plates, either MDNA-24X or MDNA-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.

Procedure

1. Prepare the reagents required for DNA library indexing.
 - a. Thaw HiFi Ultra Buffer, 5 $\times$; MM F-R Primer Mix; and MDNA-24X or MDNA-96AX index plate at room temperature.
 - b. Mix by either flicking the tube or vortexing the index plate, and then centrifuge briefly. Centrifuge the plate 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 DNA indexing

Component	Volume/reaction (µL)
Sample aliquot (from "Protocol: DNA Ligation")	6.8
HiFi Ultra Buffer, 5x	8
MM F-R Primer Mix	1.6
Nuclease-Free Water	17.1
Well from MDNA-24X or MDNA-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.

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 DNA indexing

Step	Time	Temperature (°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 DNA indexing, based on original sample input

DNA Input	FFPE sample (cycles)
40 ng	13
100 ng	12
250 ng	11

- 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 tubes/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 and Quantification". 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 starting material is a 6.8 µL sample aliquot from “Protocol: DNA Ligation”.
- 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 sample pre-treatment.
Note: Just before use, remove Buffer GE2 gDNA elimination buffer from the freezer and place on ice. After use, immediately return the tube to the freezer.
2. Prepare the sample pretreatment mix according to Table 11. 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 11. Reaction mix for sample pretreatment

Component	Volume/reaction (µL)
Sample aliquot (from “Protocol: DNA Ligation”)	6.8
Buffer GE2 gDNA elimination buffer	1.6
Total	8.4

3. Incubate the reactions in a thermal cycler according to Table 12. Use the instrument's heated lid.

Table 12. Incubation conditions for sample pretreatment

Step	Incubation temperature (°C)	Incubation time (min)
1	42	5
2	75	10
3	4	Hold

4. 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.

5. To the treated sample, prepare the reverse transcription mix according to Table 13. 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 13. Reaction mix for reverse transcription

Component	Volume/reaction (µL)
Treated Sample	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

6. Incubate the reactions in a thermal cycler according to Table 14. Use the instrument’s heated lid.

Table 14. 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

7. Add 20 μ L Nuclease-Free Water and 44 μ L QIAseq Beads and mix by vortexing or by pipetting up and down several times.
8. Incubate for 5 min at room temperature.
9. 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.
10. With the beads still on the magnetic stand, add 200 μ L 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 this second wash. Remove the ethanol with a 200 μ L pipette tip first, spin down briefly, and then use a 10 μ L 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. Ethanol carryover to the next universal PCR step will affect PCR efficiency.
13. Remove the beads from the magnetic stand, and then elute the DNA from the beads by adding 42 μ L Nuclease-Free Water.
14. Return the tubes/plate wells to the magnetic rack until the solution has cleared.
15. Transfer 40 μ L of the eluate to clean tubes/plate wells.
16. Add 44 μ L QIAseq Beads and mix by vortexing or by pipetting up and down several times.

17. Incubate for 5 min at room temperature.
18. 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.

19. With the beads still on the magnetic stand, add 200 μ L 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.
20. Repeat the ethanol wash.

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

21. 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.

22. Remove the beads from the magnetic stand, and then elute the DNA from the beads by adding 23 μ L Nuclease-Free Water.
23. Return the tubes/plate wells to the magnetic rack until the solution has cleared.
24. Transfer 20 μ L of the eluate to clean tubes/plate wells.
25. 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 starting material is the 20 μ L sample from “Protocol: RNA Reverse Transcription”.
- 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, 5 $\times$; 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 15. 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 15. Reaction mix for RNA indexing

Component	Volume/reaction (µL)
Sample aliquot (from "Protocol: RNA Reverse Transcription")	20
HiFi Ultra Buffer, 5×	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.

3. Program a thermal cycler as described in Table 16, using cycle numbers described in Table 17.

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

Table 16. Cycling conditions for RNA indexing

Step	Time	Temperature (°C)
Hold	2 min	98
2-step cycling		
Denaturation	20 s	98
Annealing	1 min	60
Extension	1 min	72
Cycle number	See Table 17	
Hold	3 min	72
Hold	∞	4

Table 17. Cycle number recommendations for RNA indexing, based on original sample input

RNA Input (ng)	FFPE sample (cycle)
80	23
200	21
500	20

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

Protocol: Target Enrichment from Indexed Whole-genome Libraries Using QIAseq xHYB Human Hybrid Capture Panels

Pool libraries and dry down the indexed library pool

Pooling of amplified, indexed whole genome libraries enables processing multiple libraries in a single hybridization-capture reaction. We recommend pooling of up to 8 indexed libraries for one hybridization capture reaction. Alternatively, hybrid-capture may be performed for a single library. The amount of indexed libraries to be pooled depends on the number of samples per pool. Make sure to only use PCR-amplified libraries for hybridization capture.

Things to do before starting

- Quantify the amplified, indexed whole genome libraries (see guidelines in “Important Notes”).
- Thaw the xHYB Probe Set, All-4-One Blocking Oligos for DNA libraries, or QIAseq N Blocking oligos for RNA libraries, and All-4-One Blocking Solution on ice. Then, pulse vortex and pulse-spin.
- To immediately proceed to hybridization capture after library pool dry-down, equilibrate the Vapor-Lock reagent to room temperature and heat the Fast Hybridization Solution to 65°C before pool evaporation is complete. Alternatively, the dried library pool can be stored at –30°C to –15°C for up to 7 days.

Procedure

- 1. For each capture pool, determine the per-library amount (ng) according to the desired pool size referring to the guidelines in Table 18 or follow the recommendations in Table 19. Make sure to not exceed 3200 ng total input per capture pool.

Table 18. Guidelines for pooling indexed libraries for hybridization capture

Library pooling guidelines

Recommended number of libraries per capture pool	8
Maximum total input per capture pool	3200 ng
Minimal input per library	200 ng

Table 19. Recommended library pooling strategy for hybridization capture

Number of indexed samples per pool	Amount of each indexed library per pool (ng)
1	200–500
2	200–500
3	200–500
4	200–500
6	200–500
8	200–400

* Do not exceed a total DNA input of 3200 ng per pool.

Note: To maintain equal representation of libraries during sequencing, use the same input amount for all libraries. Using less than 200 ng input per library may result in reduced complexity.

- 2. Calculate the volume (µL) of each amplified library needed for pooling (e.g., for a library concentration of 25 ng/µL, 16 µL are needed to add 400 ng of indexed library to the capture pool).

- Transfer the calculated volumes for each library to a 0.2 mL PCR tube/strip/plate. If the total library pool volume exceeds 180 μL , use a 1.5 mL LoBind tube. Use multiple tubes/wells if performing multiple hybridization reactions at once.

Recommendation: If the volume allows, pool libraries in a 0.2 mL PCR tube/strip or plate.

- Add the xHYB Probe Set and blocking reagents to each library pool according to Table 20.

Table 20. Complement the library pool with probes and blocking reagents

Component	Volume (μL)
Pool of indexed libraries (from step 3)	Variable
xHYB Probe Set (CGP DNA or RNA panel)*	4
One-4-All Blocking Oligos for DNA libraries or QIAseq N Blocking oligos for RNA libraries	8
One-4-All Blocking Solution	5

*Use the CGP DNA probe set for DNA library pools and the CGP RNA panel for RNA library pools.

- Completely evaporate all liquid content of the complemented library pool by using a SpeedVac system (or a similar evaporator device) equipped with an appropriate rotor for tube and/or 96-well plates. If needed, accelerate the evaporation of larger volumes by setting the temperature to 60°C.

Alternative: Evaporate all liquid content by placing the opened tube/plate in a heating block or thermal cycler with the lid opened. Evaporate at 60°C over night or until all liquid has evaporated.

- When evaporation of the complemented library pool is complete, store the dried library pool on ice and proceed immediately to step 7 (Hybridization capture protocol). Alternatively, the dried library pool can be stored at -30°C to -15°C for up to 7 days.

Hybridization capture

During the hybridization step, QIAseq xHYB Probes bind specifically to targeted library fragments present in the pool. The flexible hybridization protocol allows adjusting the duration of hybridization from as little as 30 min to 4 h. Incubating for 1 h or longer can improve the capture performance. Incubating for longer than 4 h or overnight will not enhance performance significantly, but may be considered for highest flexibility in experiment planning.

Things to do before starting

- Thaw the Vapor-Lock reagent and equilibrate to room temperature.
- Heat the Fast Hybridization Solution to 65°C in a heating block for at least 10 min and make sure all precipitate is dissolved.
- When performing the hybridization capture reaction for 1 h or less, start preparing reagents needed for binding the hybridized targets to Streptavidin Binding Beads as outlined at the beginning of the next chapter.

Procedure

7. Thaw and equilibrate the Vapor-Lock reagent to room temperature. Heat the Fast Hybridization Solution to 65°C for 10 min in a heating block. Vortex and make sure all precipitate is dissolved. Keep the Fast Hybridization Solution at 65°C until used.
8. Program a thermal cycler with a heated lid according to Table 21.

Important: The heated lid must be set to 85°C to prevent evaporation.

Table 21. Conditions for hybridization capture

Step	Incubation temperature (°C)	Incubation time
1	95	1 min
2	95	5 min
3	60	30 min–4 h*
4	60	Hold†

* Hybridization times of less than 30 min may result in reduced target region coverage. Hybridization time may be extended to overnight incubation without negative effects on specificity or uniformity.

† Program the thermal cycler to maintain 60°C after hybridization is complete. Do not allow the hybridization reaction to cool down as this negatively affects off-target rate.

9. Start the program. When the thermal cycler block reaches 95°C, pause the program in step 1.
10. Vortex the 65°C Fast Hybridization Solution and immediately transfer 20 µL to the tube/well containing the dried complemented library pool from step 6.
- Important:** Fast Hybridization Solution is viscous. Pipet slowly to ensure accurate volume transfer. For optimal volume transfer, set the pipette to 20 µL; then, slowly aspirate and dispense back into the reagent tube twice, and with the same pipette tip, slowly aspirate and transfer the entire volume to the reaction tube.
11. If the dried library pool was in a 0.2 mL PCR tube/plate, proceed to step 15.
- If the dried library pool was in a 1.5 mL LoBind tube, continue with step 12.
12. Vortex the tube for 5 s, and then incubate for 2 min at 65°C.
13. Repeat step 12 for 2 more times, and then pulse-spin the tube.

Note: A white precipitate may form. This does not affect the hybridization procedure.

14. Transfer the complete solution to a 0.2 mL PCR tube/plate and pulse-spin. Ensure that there are no bubbles present.
15. Add 30 μ L of Vapor-Lock reagent. Pulse-spin the tube to ensure all liquid is at the bottom of the tube.

Note: Vapor-Lock will form a phase on top of the hybridization reaction after pulse-spin. This does not affect the hybridization procedure.

16. Transfer the tube/plate to the thermal cycler and continue the program at 95°C.

Important: Make sure the heated lid is set to 85°C and the tube is tightly sealed with fresh lids/foil. Make sure to use the correct lid spacers for your thermal cycler.

17. Leave the hybridization reaction in the thermal cycler until ready for binding of captured library fragments to Streptavidin beads. Continue with the protocol “Bind hybridized targets to Streptavidin beads” at least 30 min before hybridization is complete.

Alternatively, the hybridization reaction may remain in the thermal cycler over night to continue with binding the hybridized targets to Streptavidin beads the next day.

Important: Do not allow the hybridization reaction to cool to less than 60°C after the program is complete.

Bind hybridized targets to Streptavidin beads

In this step, the biotin-coupled xHYB Probes and their hybridized targets will be captured on Streptavidin beads. Nonspecific targets will be washed away during a series of washing steps. Start preparing the Streptavidin Binding Beads and wash buffers at least 30 min before hybridization is complete.

Things to do before starting

- Preheat a heating block for 1.5 mL tubes to 70°C.
- Preheat a heating block for 1.5 mL tubes to 48°C.
- Heat the whole reagent bottles of the Wash Buffer A and Wash Buffer B to 48°C and mix thoroughly. Leave both reagent bottles at 48°C until needed.
- Heat the whole reagent bottle of the Post Capture Binding Buffer to 48°C and mix thoroughly. Then leave the reagent bottle at room temperature until needed.
- Equilibrate Streptavidin Binding Beads to room temperature for at least 30 min.

Procedure

18. Per reaction, aliquot 450 µL Wash Buffer A and preheat to 70°C.

19. Per reaction, aliquot 700 µL Wash Buffer B and preheat to 48°C.

Important: For steps 18 and 19, do not aliquot chilled buffers. Heat the entire reagent bottles of Wash Buffer A and Wash Buffer B to 48°C and mix thoroughly. Then, prepare aliquots.

Note: Leave the wash buffer aliquots in the heating blocks until needed.

20. Homogenize the pre-equilibrated Streptavidin Binding Beads by vortexing.

21. Transfer 100 µL homogenized Streptavidin Binding Beads to a new 1.5 mL LoBind tube. Prepare 1 tube for each hybridization reaction.

22. Add 200 µL Post Capture Binding Buffer and mix by pipetting.

Important: Do not use chilled Post Capture Binding Buffer. Make sure to heat the buffer to 48°C before use and mix well. Then allow the buffer to cool to room temperature.

23. Incubate the beads on a magnetic stand for 1 min or until the solution is clear, and then remove and discard the clear supernatant without disturbing the pellet. Remove from the magnetic stand.
24. Repeat steps 22 and 23 for 2 more times for a total of 3 washing steps.
25. After removing the supernatant from the third wash, add a final 200 μ L Post Capture Binding Buffer and homogenize the beads by vortexing. Keep the homogenized bead slurry at room temperature until needed.
26. After the hybridization reaction is complete, open the lid of the thermal cycler, open the tube, and swiftly transfer the complete volume including Vapor-Lock into the tube containing the bead slurry from step 25. Mix well by pipetting.

Important: Do not allow the hybridization reaction to cool to less than 60°C before transferring to the Streptavidin Binding Beads. Rapid transfer directly from the thermal cycler is essential for minimizing the off-target rate.

27. Incubate the mixture at room temperature for 30 min on a shaker, rotator, or similar device to prevent the beads from settling.

Alternatively, incubate the mixture at room temperature for 30 min, while agitating in 5-min intervals to prevent beads from settling.

Note: Aggressive mixing is not required. Do not vortex.

28. Pulse-spin the tube to ensure all liquid is at the bottom of the tube.
29. Incubate the beads on a magnetic stand for 1 min or until the solution is clear, and then remove and discard the clear supernatant including Vapor-Lock without disturbing the pellet.

Note: Small amounts of Vapor-Lock may remain after removal of supernatant and throughout each wash step. This does not affect the final capture product.

30. Remove the tube from the magnetic stand, and add 200 μ L of preheated 70°C Wash Buffer A. Mix by pipetting.
31. Incubate for 5 min at 70°C.
32. Place the tube on a magnetic stand for 1 min or until the solution is clear, and then remove and discard the clear supernatant without disturbing the pellet.

Remove the tube from the magnetic stand, and add another 200 μ L of preheated 70°C Wash Buffer A. Mix by pipetting.
33. Incubate for 5 min at 70°C.
34. Pulse-spin the tube and transfer the entire volume, including the beads to a new 1.5 mL LoBind tube.
35. Place the tube on a magnetic stand for 1 min or until the solution is clear, and then remove and discard the clear supernatant without disturbing the pellet.
36. Remove the tube from the magnetic stand, and add 200 μ L of preheated 48°C Wash Buffer B. Mix by pipetting.
37. Pulse-spin the tube, and then incubate for 5 min at 48°C.
38. Place the tube on a magnetic stand for 1 min or until the solution is clear, and then remove and discard the clear supernatant without disturbing the pellet.
39. Repeat steps 36 to 38 for 2 more times for a total of 3 washes.
40. After the final wash, use a 10 μ L pipette to remove all residual supernatant. Proceed immediately to the next step.

Important: Do not allow the beads to dry.
41. Remove the tube from the magnetic stand, and add 45 μ L RNase-free water. Homogenize by pipetting, and store the bead slurry on ice.
42. Proceed directly to “Postcapture amplification”.

Alternatively, aliquot the bead slurry into two 22.5 µL aliquots and store at –30°C to –15°C for later use.

Important: Do not discard the beads. The capture product is still bound to the Streptavidin beads.

Postcapture amplification

Captured targets bound to Streptavidin Binding Beads will be amplified using QIAGENs Post Hybrid-Capture PCR Mix. The amplified hybrid capture library will be purified and is ready for QC and sequencing on Illumina instruments.

Things to do before starting

- Prepare fresh 80% ethanol.
- If stored at –30°C to –15°C, thaw a 22.5 µL aliquot of the Streptavidin Binding Bead slurry on ice (from step 42).
- Equilibrate Agencourt AMPure XP beads to room temperature for 20–30 min before use.
- Thaw Post Hybrid-Capture PCR Mix and Primer Mix Illumina Library Amplification on ice. Once reagents are thawed, mix by pulse-vortexing to avoid any localized concentrations. Briefly spin down vortexed reagents before use.

Procedure

43. Program a thermal cycler with a heated lid according to Table 22 and Table 23.

Table 22. Postcapture amplification conditions

Time	Temperature (°C)	Number of cycles
2 min	98	1
20 s	98	7–13 cycles*
30 s	60	
30 s	72	
1 min	72	1
∞	4	Hold

* See Table 23 for the optimal cycle numbers depending on input amount and xHYB panel used.

Table 23. Number of postcapture amplification cycles

xHYB Panel	<500 ng capture input	500 ng – 2.0 µg capture input	>2.0 µg capture input
CGP DNA Panel	12 cycles	11 cycles	10 cycles
CGP RNA Panel	13 cycles	12 cycles	11 cycles

44. Prepare the reaction mix on ice according to Table 24. Mix the components in a PCR tube or 96-well PCR plate.

Table 24. Reaction mix for postcapture amplification

Component	Volume (µL)
Bead slurry with captured targets (from step 41)	22.5
Post Hybrid-Capture PCR Mix	25
Primer Mix Illumina Library Amplification	2.5
Total reaction volume	50

45. Transfer the PCR tube or plate to the thermal cycler and start the program.
46. Once PCR is complete, add 75 μL of resuspended QIAseq beads to each reaction (50 μL) and pipet up and down thoroughly to mix.
47. Incubate the mixture for 5 min at room temperature. Pellet the beads on a magnetic stand, and carefully discard the supernatant.
48. Wash the beads by adding 200 μL of 80% ethanol. Pellet the beads on the magnetic stand and discard the supernatant. Repeat the wash once for a total of 2 ethanol washes. Remove as much excess ethanol as possible.
49. Incubate on the magnetic stand for 5–10 min or until the beads are dry. Overdrying of beads may result in lower DNA recovery. Remove from the magnetic stand.
50. Elute by resuspending in 32 μL of Buffer EB, 10 mM Tris-Cl, pH 8.0, or RNase-free water. Incubate at room temperature for 2 min.
51. Pellet the beads on the magnetic stand. Carefully transfer 30 μL of the clear supernatant into a new tube.
52. Assess the quality and size distribution of the library using a capillary electrophoresis device such as QIAGEN QIAxcel Advanced or Agilent Bioanalyzer 2100. Check for the expected size distribution (see Figure 2). For best sequencing results, quantitate the captured library using a qPCR assay such as the QIAseq Library Quant Assay Kit or compatible method.

Note: The library should show a distribution similar to the whole genome libraries used for the respective capture pool.
53. The purified library can be safely stored at -30°C to -15°C in a DNA LoBind tube until ready to use for sequencing.

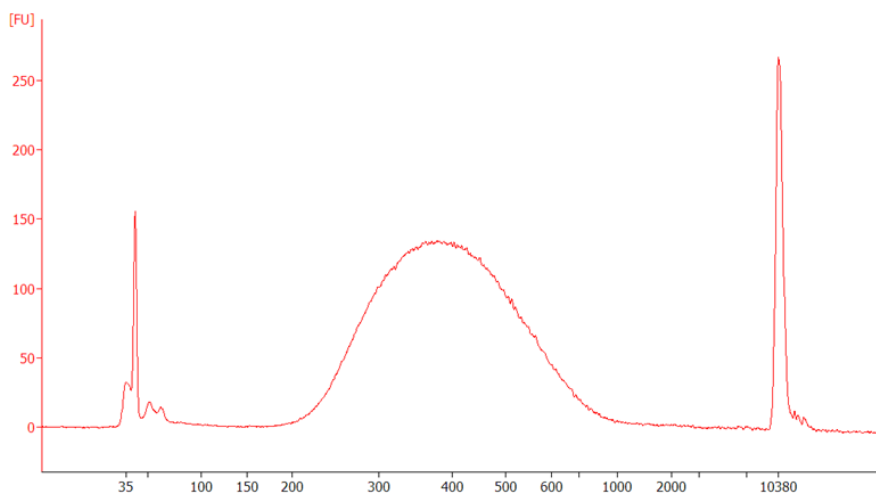


Figure 2. Electrophoresis trace data of a final hybrid capture library pool.

Recommendations: Library QC and Quantification

NGS Library QC

Perform QC with the QIAxcel Connect or similar instrument. Check for the correct size distribution of library fragments (~400–500 bp median size) and for the absence of adapters or adapter–dimers (~<200 bp) (Figure 3).

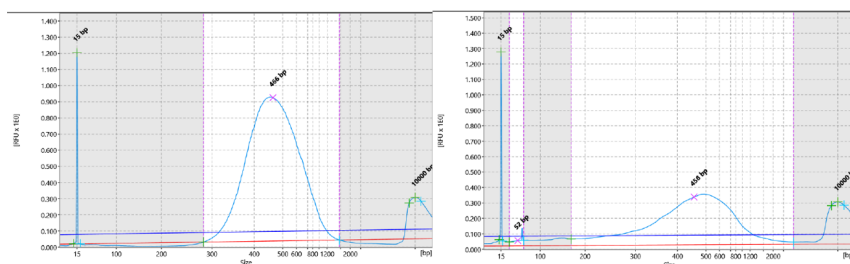


Figure 3. QIAseq Multimodal WGS DNA (left) and WTS RNA (right) Libraries.

Preferred library quantification method

Quantify the library using Nanodrop® or QIAxpert® if proceeding to “Hybrid Capture”, or qPCR-based method if proceeding to sequencing.

Sequencing Guidelines

Pooling and multiplexing guidelines

Depending on the sequencing instrument used, combining multiple hybridization capture pools may be advisable for cost efficient sequencing. Before combining equal volumes of hybrid capture pools, quantify each pool and dilute to a defined concentration (e.g., 5 nM). To determine the total number of samples to be sequenced in a single sequencing run, consider specifications of the sequencing instrument, the safe target cluster density, target region size of the used xHYB panel and the desired coverage per sample. Please refer to Table 25a and Table 25b for initial multiplexing guidelines per sequencing run to achieve a minimal average target region coverage of 1200x for DNA Library. Multiplexing and loading may need to be adjusted based on your local conditions like the quantification method used and your preferences for target cluster density.

xHYB panels achieve a very high coverage uniformity throughout the target region with a typical Fold-80 base penalty in the range of 1.3–1.5 or better. This means that the average target region coverage needs to be increased by approximately 1.4 to raise 80% of non-zero-coverage target bases to the mean coverage level. When planning the sequencing pool, consider approximately 1.3- to 1.5-fold of the desired minimal coverage to reliably obtain that coverage throughout the target region.

Table 25a. Multiplexing guidelines for Illumina instruments using 2 × 149 bp paired-end sequencing to achieve a minimal average target coverage of 1200x for DNA Library. Actual coverage may vary depending on input DNA quality and the sequencing instrument used

	# (DNA library+RNA library)	# (DNA library)	# (RNA library)*
DNA library reads (M)	75	75	–
Reads for RNA library (M)	50	–	50

*When sequencing an RNA library alone (without DNA), we recommend adding 20% pHIX.

Table 25b. Multiplexing guidelines for Illumina instruments using 2 × 149 bp paired-end sequencing to achieve a minimal average target coverage of 1200× for DNA Library. Actual coverage may vary depending on input DNA quality and the sequencing instrument used

Platform and Sequencing Chemistry

Instrument	Version	Read length	Capacity (paired-ends reads)	Capacity (M)	# (DNA library+RNA library)	# (DNA library)	# (RNA library)*
NextSeq® 500/550	Mid output	2 × 149 bp	260 M	260	2	3	5
NextSeq 500/550	High output	2 × 149 bp	800 M	800	6	10	16
NextSeq 1000/2000	P1	2 × 149 bp	200 M	200	1	2	4
NextSeq 1000/2000	P2	2 × 149 bp	800 M	800	6	10	16
NextSeq 1000/2000	P3	2 × 149 bp	2400 M	2400	19	32	48
NovaSeq™ 6000	SP (2 lanes per flow cell)	2 × 149 bp	1.6 B	1600	12	21	32
NovaSeq 6000	S1 (2 lanes per flow cell)	2 × 149 bp	3.2 B	3200	25	42	64
NovaSeq 6000	S2 (2 lanes per flow cell)	2 × 149 bp	8.2 B	8200	65	109	164
NovaSeq 6000	S4 (4 lanes per flow cell)	2 × 149 bp	20 B	20,000	160	266	400
NovaSeq X Series	1.5 B	2 × 149 bp	3.2 B	3200	25	42	64
NovaSeq X Series	10 B	2 × 149 bp	20 B	20,000	160	266	400
NovaSeq X Series	25 B	2 × 149 bp	52 B	52,000	416	693	1040

Table 25b. Multiplexing guidelines for Illumina instruments using 2 × 149 bp paired-end sequencing to achieve a minimal average target coverage of 1200× for DNA Library. Actual coverage may vary depending on input DNA quality and the sequencing instrument used (continued)

Platform and Sequencing Chemistry

Instrument	Version	Read length	Capacity (paired-ends reads)	Capacity (M)	# (DNA library+RNA library)	# (DNA library)	# (RNA library)*
AVITI/AVITI LT	Low Output (per flow cell)	2 × 149 bp	500 M	500	4	6	10
AVITI/AVITI LT	Midium Output (per flow cell)	2 × 149 bp	1 B	1000	8	13	20
AVITI	High Output (per flow cell)	2 × 149 bp	2 B	2000	16	26	40

*When sequencing an RNA library alone (without DNA), we recommend adding 20% pHix.

Sequencing

- Always ensure that libraries have been quantified using QIAseq Library Quant Assay or a compatible method to enable equal library representation within the sequencing pool and exact pool concentrations for optimal flow cell loading and best sequencing performance.
- For complete instructions on how to denature sequencing libraries and set up a sequencing run, please refer to the system-specific Illumina documents.
- Depending on the adapter type used for library preparation, please refer to the respective QIAseq Library handbook for detailed information on QIAseq Combinatorial and Unique Dual-Index Adapters.
- Editable and ready to use sample sheets, including all sample indexes of QIAseq Multimodal DNA/RNA UDI are available for download at www.qiagen.com

- A description of run setup for Illumina instruments and definition files for QIAseq Multimodal DNA/RNA UDI to be used for Illumina Local Run Manager are available at www.qiagen.com
- Sequencing on the NextSeq, HiSeq X®, HiSeq® 3000/4000 systems follow a different dual-indexing workflow than other Illumina systems. If manually creating sample sheets on these instruments, enter the reverse complement of the i5 index adapter sequence. If using BaseSpace®, or Local Run Manager to plan the run, the software will automatically reverse complement index sequences when necessary.

Sequencing setup on Illumina instruments

The following guidelines outline the most important settings for Illumina instruments. More detailed instructions on how to configure a run and how to create a sample sheet can be found in the “Product Resources” at www.qiagen.com

- Read Type: **Paired End**
- Index Reads: **2**
- Enable Adapter Trimming
- Cycles:
 - Read 1 and Read 2: **149**
 - Index 1 and Index 2: **10**

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 at our Technical Support Center: www.qiagen.com/FAQ/FAQList.aspx (for contact information, visit www.qiagen.com).

Comments and suggestions	
Low CGP hybrid capture library yields	
a. Suboptimal hybrid capture input quality or amount	Make sure to use high-quality indexed whole genome libraries as input that have been quantified using a fluorometric method or capillary electrophoresis.
b. Hybridization time too short	Hybrid capture input amounts smaller than 1000 ng may result in decreased library yields when using hybridization times of 30 min or less. Increase hybrid capture input, hybridize for 1 h or longer.
c. Inappropriate amount of Streptavidin Binding Beads	Insufficient Exome library yield may be a result of deficient postcapture amplification. Streptavidin Binding beads may inhibit postcapture amplification. Ensure to use no more than 100 µL of Streptavidin Binding beads per hybridization capture reaction. Only amplify half of the final bead slurry per PCR.
d. Inappropriate hybridization or washing temperature	Hybridization capture performs optimally at a very narrow temperature range. Ensure your thermal cycler and heating blocks are calibrated to accurately reach the intended temperature for the hybridization reaction and washing steps.
Decreased target region specificity or uniformity of CGP DNA libraries	
a. Incorrect use of Blocking Oligos or Blocking Solution	Only use QIAseq One-4-All Blocking Oligos and One-4-All Blocking Solution as outlined in the protocol.
b. Wrong library adapter type used	One-4-All Blockers delivered with the kit are designed for Illumina TruSeq compatible adapters, but not for other adapter types. N-Blockers are designed for Illumina Nextera compatible adapters.
c. Suboptimal hybridization and washing temperature	Check thermal cycler programs and make sure thermal cyclers and heating blocks are properly calibrated. For some heating blocks, it may be necessary to adjust the washing temperatures by 1°C or 2°C to match the optimal washing conditions.

Comments and suggestions

- | | |
|--|--|
| d. Suboptimal hybridization and washing conditions | Follow the protocol for hybridization and washing steps. Take care not to confuse the usage order of Wash Buffers A and B. Ascertain that Fast Hybridization Solution is heated to 65°C before use and neither the Fast Hybridization Solution nor the Wash Buffers show precipitates when aliquoting. |
|--|--|

Unequal read representation of libraries within a hybrid capture pool

- | | |
|---|--|
| a. Unequal hybrid capture input | Use equal amounts of indexed whole genome libraries to be pooled for a hybrid capture reaction. Quantify indexed whole genome libraries before hybrid capture by using a capillary electrophoresis method such as the QIAxcel Advanced System or the Agilent 2100 Bioanalyzer using High Sensitivity DNA chips. |
| b. Suboptimal library enrichment conditions | Do not use PCR-free whole genome libraries. Always follow the library enrichment protocol outlined in the respective library preparation protocol using QIAGEN HiFi PCR Master Mix for library enrichment. Make sure input libraries are not overamplified. Quantify indexed whole genome libraries using the QIAGEN Library Quant Assay and assess library quality via capillary electrophoresis. |

Ordering Information

Product	Contents	Cat. no.
QIAseq xHYB CGP DNA Panel (24)	xHYB probe panel for DNA target enrichment for comprehensive genomic profiling, fixed panel for 24 samples	333122
QIAseq xHYB CGP DNA Panel (96)	xHYB probe panel for DNA target enrichment for comprehensive genomic profiling, fixed panel for 96 samples	333125
QIAseq xHYB CGP RNA Panel (24)	xHYB probe panel for RNA target enrichment for comprehensive genomic profiling, fixed panel for 24 samples	334122
QIAseq xHYB CGP RNA Panel (96)	xHYB probe panel for RNA target enrichment for comprehensive genomic profiling, fixed panel for 96 samples	334125
QIAseq xHYB CGP DNA/RNA Panels (24)	xHYB probe panels for DNA and RNA target enrichments for comprehensive genomic profiling, fixed panel for 24 samples	335122
QIAseq xHYB CGP DNA/RNA Panels (96)	xHYB probe panels for DNA and RNA target enrichments for comprehensive genomic profiling, fixed panel for 96 samples	335125
QIAseq xHYB Reagent Kit (24)	Reagents required for hybrid capture-based target enrichment. The kit contains 3 hybrid capture reactions, which are typically sufficient for 24 samples.	333430
QIAseq xHYB Reagent Kit (96)	Reagents required for hybrid capture-based target enrichment. The kit contains 12 hybrid capture reactions, which are typically sufficient for 96 samples.	333195

QIAseq library kits for the generation of indexed whole genome and whole transcriptome libraries for use in hybridization capture

QIAseq Multimodal DNA/RNA Lib Kit (24)	Reagents (except indexes) for multimodal (DNA and RNA) library preparation of 24 samples: 24 DNA and 24 RNA libraries	334842
QIAseq Multimodal DNA/RNA Lib Kit (96)	Reagents (except indexes) for multimodal (DNA and RNA) library preparation of 96 samples: 96 DNA and 96 RNA libraries	334845
QIAseq Multimodal DNA/RNA UDI (24)	DNA and RNA indexes for 24 samples for a total of 48 reactions	334852
QIAseq Multimodal DNA/RNA UDI (96)	DNA and RNA indexes for 96 samples for a total of 192 reactions	334855

Product	Contents	Cat. no.
QIAseq N Blocking Oligos (12)	Oligos designed to bind and block all Illumina Nextera-compatible NGS adapters for 12 reactions of hybridization capture	334871
QIAseq Library Quantification Kits for use with Illumina instruments		
QIAseq Library Quant Assay Kit	Laboratory-verified forward and reverse primers for 500 × 25 µL reactions (500 µL); DNA Standard (100 µL); Dilution Buffer (30 mL); (1.35 mL × 5) GeneRead® qPCR SYBR® Green Mastermix	333314

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 Support or your local distributor.

Document Revision History

Date	Description
04/2025	Initial release.
04/2026	Correction in Kit Contents. Updated "Nucleic Acid Fragmentation", "RNA Polyadenylation", "DNA Ligation", "DNA Library Indexing", "RNA Reverse Transcription", and "RNA Library Indexing".

Limited License Agreement for QIAseq® xHYB CGP DNA/RNA Panel

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®, QIAamp®, AllPrep®, FastSelect®, QCI®, PAXgene®, QIAxcel® (QIAGEN Group); Agilent® B (Agilent Technologies); LoBind® (Eppendorf AG); Nextera®, Illumina®, iSeq®, TruSeq®, HiSeq®, MiniSeq®, MiSeq®, NextSeq®, NovaSeq® (Illumina, Inc.); MagneSphere® (Promega Corporation); DynaMag™, NanoDrop®, Quant-iT™ (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.

06/2026 HB-3690-002 © 2026 QIAGEN, all rights reserved.

Notes.

Notes.

Notes.

