

Biomek i7 Hybrid Liquid Handler Automated Method for QIAseq® miRNA UDI Library Kit



2022-GBL-EN-100479-v1

Notices

Biomek Automated Workstations are not intended or validated for use in the diagnosis of disease or other conditions.

QIAseq® miRNA Library Kits are intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals. QIAseq® is a trademark of QIAGEN Group. QIAseq® miRNA Library Kits are sold by QIAGEN Group.

© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. This protocol is for demonstration only; and is not validated by Beckman Coulter.

QIAseq® miRNA UDI Library Kit

- Qiagen's QIAseq® miRNA UDI Library kit enables Sample to Insight®, precision next-generation sequencing (NGS) of mature miRNAs on Illumina® NGS instruments.
- QIAseq® miRNA UDI Library Kit does not require gel purification, excision, and elution, which substantially reduces the required hands-on time and noticeably shortens the length of the whole workflow.
- The QIAseq® miRNA UDI Library Kit uses mature miRNA and contains the following steps:
 1. 3' ligation
 2. 5' ligation
 3. cDNA synthesis
 4. cDNA cleanup
 5. library amplification
 6. library cleanup

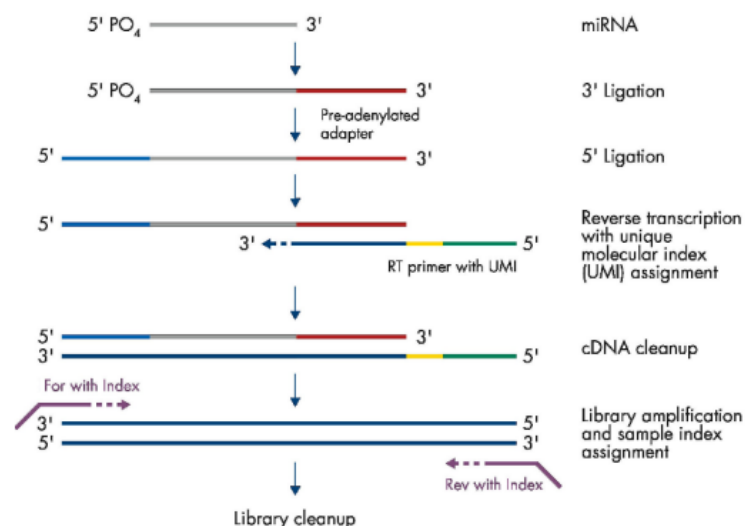


Figure 1: QIAseq® miRNA UDI Library Kit Workflow

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Automated Method Description:

- Sample number ranging from 1-96
- On-Deck or Off- Deck Thermocycler
- RNA sample inputs of 500ng, 100ng, 10ng, 1ng, or serum/plasma
- User defined starting well position for UDI index plate
- Optimization of PCR cycles



Figure 2: Biomek i7 Hybrid Liquid Handler

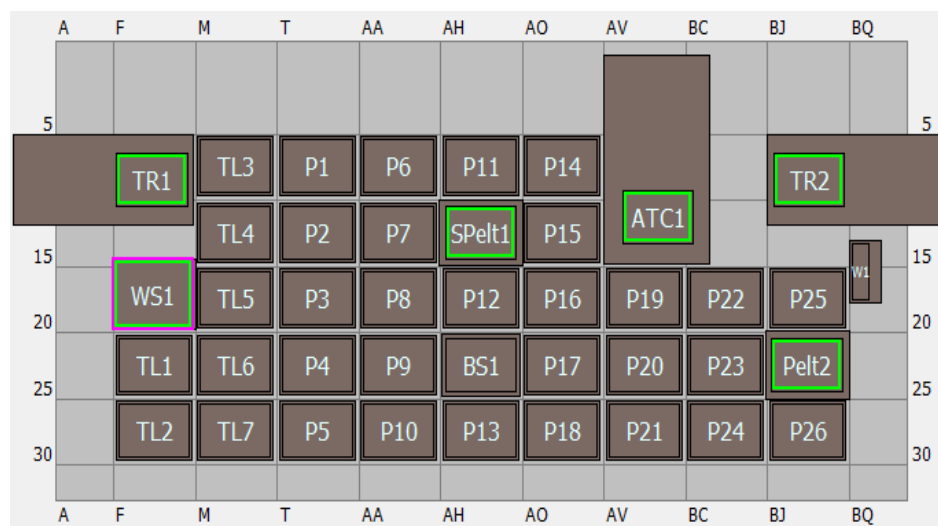


Figure 3: Biomek i7 Hybrid Liquid Handler Deck Layout

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Part Number	Qty	Manufacturer	Description
B87585	1	Beckman Coulter	Biomek i7 Hybrid (MC + Span-8) with Enclosure
	1	Beckman Coulter	Pipetting head, MC-96, 5 - 1200 µL
	1	Beckman Coulter	Pipetting head, Span-8, 1mL
C02613	1	Beckman Coulter	Mobile Workstation, i7
719366	2	Beckman Coulter	Device Controller I/O Box
B87483	2	Beckman Coulter	ALP, Trash
C02867	7	Beckman Coulter	ALP, Tip Load 1x1
B87477	23	Beckman Coulter	ALP, Static 1x1
B87478	1	Beckman Coulter	ALP, 1x3 static
A93942	1	Beckman Coulter	ALP, Shaking Peltier Kit
C05041	1	Beckman Coulter	Kit, Integration - Bioshake on i-Series
C02750	2	Beckman Coulter	Vibration isolation mounting plate
A93938	1	Beckman Coulter	ALP, Static Peltier Kit
B87485	1	Beckman Coulter	ALP, Mounting Plate
B87689	1	Beckman Coulter	Wash station, MC-96 pipette tips
719654	1	Beckman Coulter	Wash station, Span-8 pipette tips, Passive
C05021	1	Beckman Coulter	Kit, Integration, Deck - ATC on i-Series
C05753	1	Beckman Coulter	Kit, Installation, i-series Workstation (including Biomek Config Tool Software)
Contact Manufacturer	1	Thermo Fisher	Automated Thermocycler (ATC), 96 well block

Table 1: Biomek i7 Hybrid Liquid Handler Part List

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

The QIAseq® miRNA UDI Library Kit automation method for the Biomek i7 Liquid Handler can be completed as 2 sections run independently or combined for the complete workflow.

The method utilizes a single user interaction for the guided deck setup and can process up to 96 samples. The completion of the method will result in sequence-ready libraries.

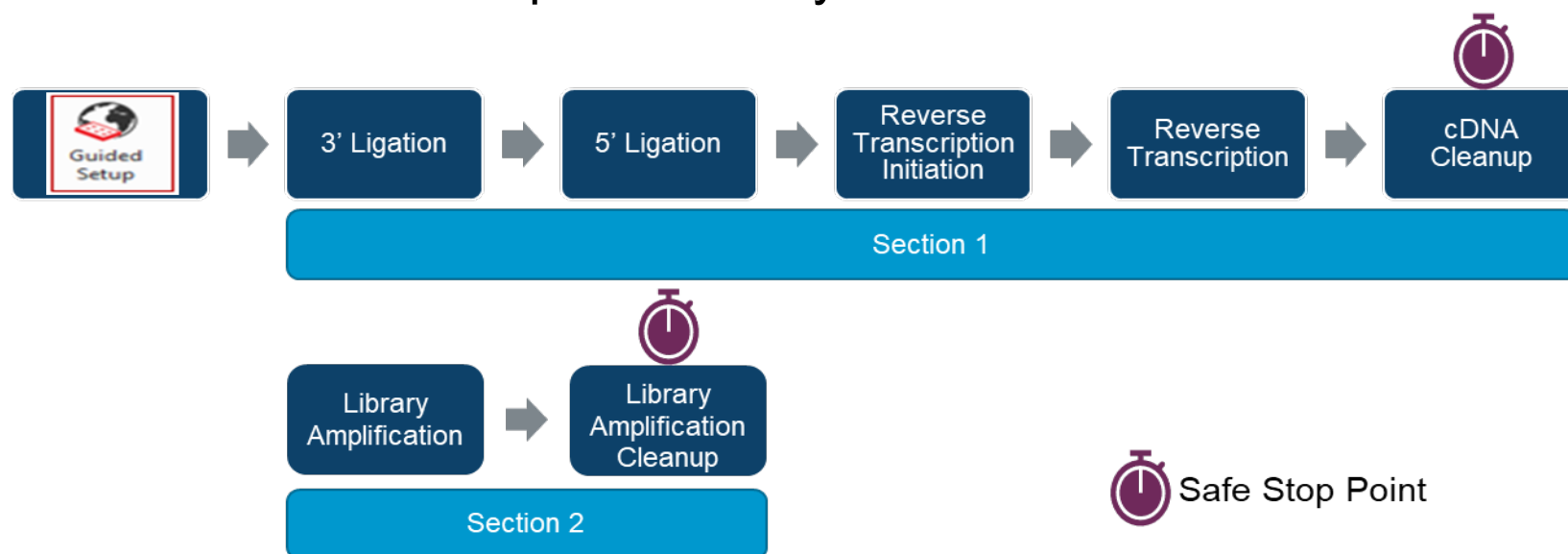


Figure 4: QIAseq® miRNA UDI Library Kit automated workflow for the Biomek i7 Hybrid Liquid Handler

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Section Information	Timing for 24 Samples	Timing for 96 Samples
Instrument Setup	10 minutes	15 minutes
Section 1: 3' Ligation, 5' Ligation, Reverse Transcriptase, and Cleanup	6 hours	7 hours
Section 2 : Library Amplification and Cleanup	2 hours, 30 minutes	3 hours
Total Run Time	8 hours, 40 minutes	10 hours, 15 minutes
*Total time estimates do not include reagent thawing and preparation. One user interaction is needed for the instrument setup.		

Table 2: QIAseq® miRNA UDI Library Kit automation timing estimates

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Demonstrated Method Interface

Biomek Method Launcher

1. Makes method available outside of the editor
2. Protects against accidental changes from multiple users

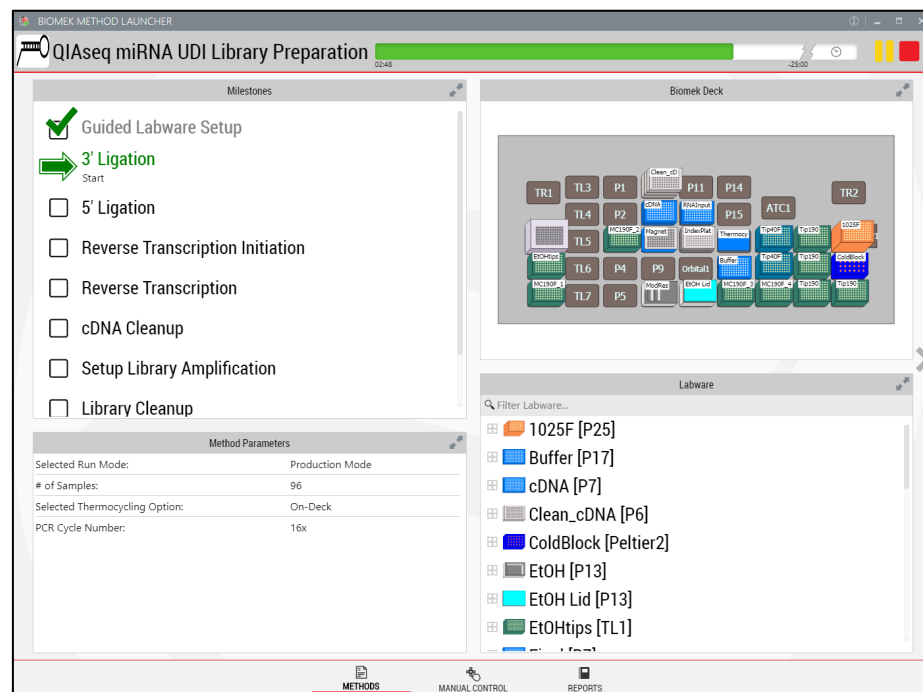


Figure 5: QIAseq® miRNA UDI Library Kit automated workflow for the Biomek i7 Hybrid Liquid Handler as shown on the Biomek Method Launcher interface

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Demonstrated Method Interface

Method Option Selector (MOS)

Dynamic, HTML-driven interface allows the user to select a number of options to configure the workflow as desired.

Full walk away capability from beginning to end of the protocol with multiple restart locations.

Beckman Coulter

QIAseq® miRNA UDI Library Kit

Optimized for Biomek i7 Dual Hybrid Automated by Beckman Coulter

Method Parameters

Number of Samples: (1-96) Samples

Sample Type:

Select Ethanol Drying Time (minutes): (10-25)

Use On-Deck Thermocycler?

Method Options

☒ 3' Ligation, 5' Ligation, Reverse Transcription, and cDNA Cleanup

☒ Library Amplification and Cleanup

Index Plate Starting Position:

Select Number of PCR Cycles:

BECKMAN COULTER

Figure 6: QIAseq® miRNA UDI Library Kit automated workflow for the Biomek i7 Hybrid Liquid Handler Method Option Selector

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Demonstrated Method Interface

MOS options that enable flexibility in configuring a run:

- 1. Number of Samples:** Allows the user to manually enter the number of samples (1-96) to be processed.
- 2. Sample Type:** Options include 500ng, 100ng, 10ng, 1ng, or serum/plasma.
- 3. Ethanol Drying Time:** User can define the drying time for Ethanol depending on lab conditions. Options range from 10-25 minutes.
- 4. Use On-Deck Thermocycler:** Results in a drop-down menu of either “yes” or “no” to match the deck layout of the user’s Biomek i7 hybrid Automated Workstation.
- 5. Index Plate Starting Position:** Drop-down menu that allows the user to specify the starting location within the 96-well Index plate.
- 6. Select Number of PCR Cycles:** Allows the user to set the number of PCR cycles run for Library Amplification. The number of cycles are supported from 12-26.

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Demonstrated Method Interface

Guided Labware Setup (GLS) and Reagent Calculation

- Provides clear instructions for setup of labware
- Provides calculated volumes of reagents based on the number of samples processed

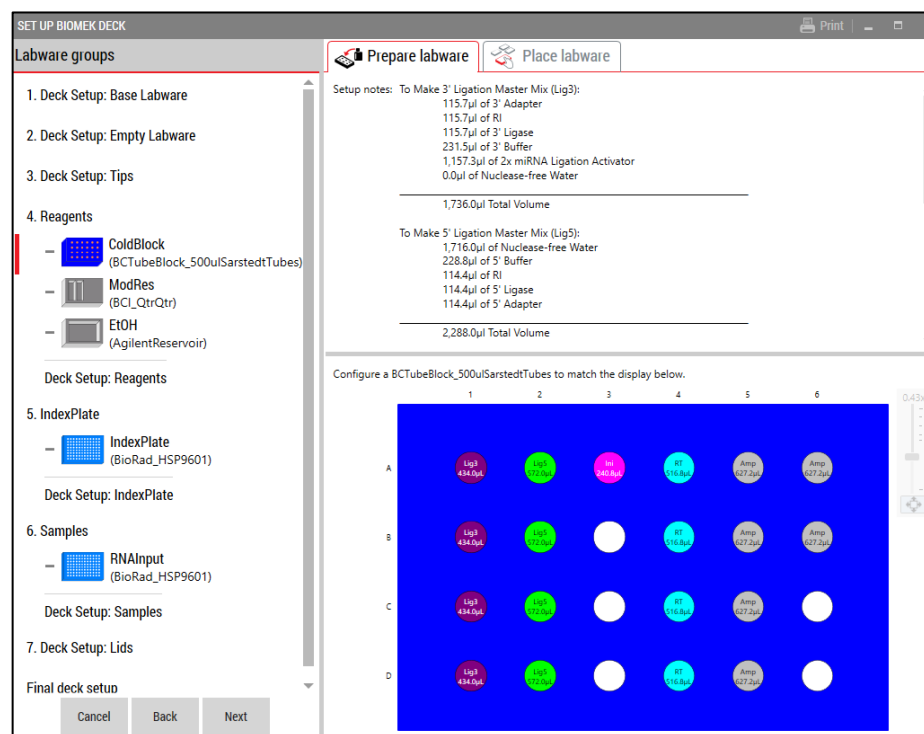


Figure 7: Guided Labware Setup for QIAseq® miRNA UDI Library Kit on the Biomek i7 Hybrid Liquid Handler

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

Automation runs were performed using 24 technical replicates of Human XpressRef Universal Total RNA (Qiagen, 338112) at 100 ng and 10 ng. Following library preparation, the libraries were analyzed using High Sensitivity D1000 ScreenTapes (Agilent, 5067-5584, 5067-5585) on the Agilent TapeStation Instrument for library size, concentration, and quality.

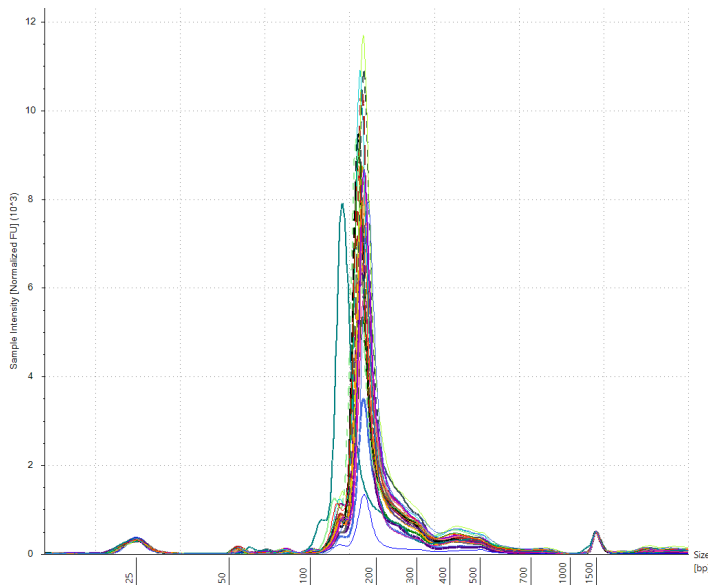


Figure 8: XpressRef 100ng libraries from QIAseq® miRNA UDI Library Kit on the Biomek i7 Hybrid Liquid Handler

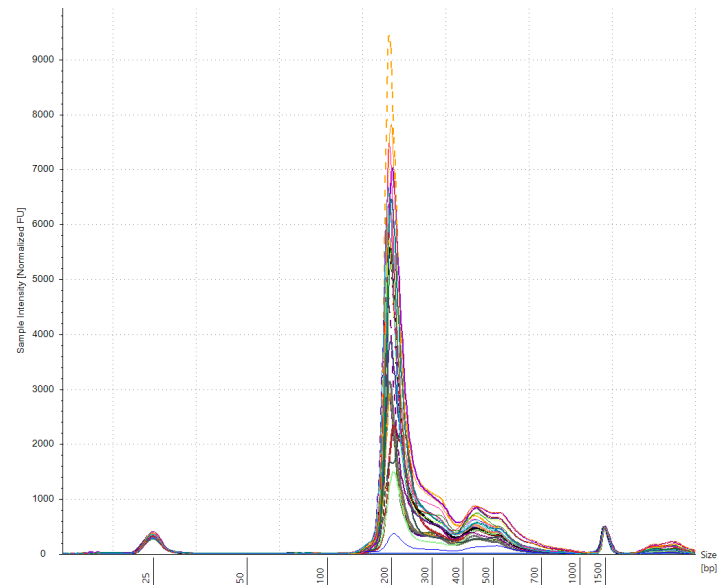


Figure 9: XpressRef 10ng libraries from QIAseq® miRNA UDI Library Kit on the Biomek i7 Hybrid Liquid Handler

Biomek i7 Hybrid Liquid Handler automated method for QIAseq® miRNA UDI Library Kit

High Sensitivity DNA analysis (Agilent TapeStation) of 24 technical replicates of Human XpressRef Universal Total RNA (Qiagen, 338112) was performed. The 100 ng libraries demonstrated an average library size of 177bp and average concentration of 6685 pg/μL. The 10 ng libraries demonstrated an average size of 201bp and average concentration of 4689 pg/μL.

Sample Type	Average Size (bp)	Average Concentration (pg/μL)
XpressRef (100 ng)	177	6685
XpressRef (10 ng)	201	4689

Table 3: Average size and concentration of 100ng and 10ng QIAseq® miRNA UDI libraries from the Biomek i7 Hybrid automated method.

Intellectual Property

This document contains confidential and proprietary information owned by Beckman Coulter, Inc.

Any disclosure, use, copying or distribution of this information without the express written consent of Beckman Coulter, Inc. is strictly prohibited.