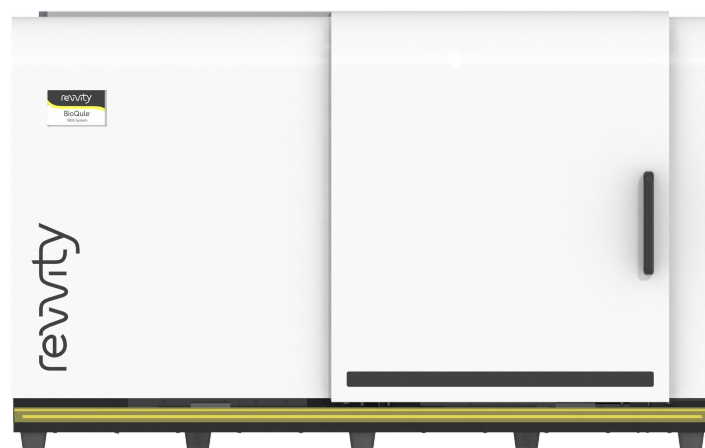




## Assay Guide

# BioQule™ Illumina® DNA Prep Method

Copyright 2022 - 2024 Revvity Inc. All rights reserved.  
PN: PR-00006 Rev E  
Publication Date: July 19, 2024



# Table of Contents

1	Introduction	3
1.1	Overview	3
1.2	Storage and Stability	3
1.3	Product Use	4
1.4	Warnings and Precautions	4
1.5	Prior to Starting	4
2	Contents	5
2.1	Kit Contents	5
2.2	Additional Equipment, Reagents and Labware	5
3	Planning the Run	6
3.1	Workflow and Time Required	6
3.2	Input DNA Requirements	6
3.3	Sequencing Recommendations and Guidelines	7
3.4	Data Analysis	7
3.5	Library Storage	7
4	Procedure	8
4.1	Reagent Plate Setup	8
4.2	BioQule Run Setup	11
5	Troubleshooting	14
5.1	Low Volume of Library Generated	14
5.2	Low Yield Library Generated	15
5.3	High Adapter Dimer	15
6	Appendix	16
6.1	Index (UDI) Sequences	16
7	Technical Assistance	21
8	Revision History	21

# 1 Introduction

## 1.1 Overview

The BioQule™ NGS system is an innovative automated platform that simplifies most commercially available Next Generation Sequencing workflows with the push of a button. It employs simplified micro and macroscale geometries to efficiently perform DNA extraction, library preparation, and library quantification. DNA libraries prepared using the platform meet high-quality standards regarding coverage bias, yield, and fragment size. It is an open system that can integrate and automate different NGS workflows, including Illumina, Element Biosciences, and Oxford Nanopore Technologies NGS workflows.

BioQule is ideal for laboratory setups where high-throughput robotic liquid handlers are not feasible. It has been designed so that no automation experience is required to prepare NGS libraries. A user only has to prepare the pre-plated reagent plate, place the cartridge and plate into the instrument, close the door, and start the instrument. Hence, the BioQule offers full walkaway automation (80% reduction in hands-on time) and is easy to use (only 3 touch points). This avoids day-to-day and sample-to-sample human error (>99% pass rate).

### Features:

- Low input requirement for genomic DNA down to 40 ng/sample.
- Complete library prep solution, including size selection beads and fluorescence measurements for quantifying libraries.
- Robust genome coverage and reliable performance with sequencing bias mitigation.
- Functionally tested with Illumina sequencing platform.

### BioQule™ Illumina® DNA Prep Specifications:

Input Type:	gDNA
Input Amount:	40 ng - 500 ng
Number of Reactions:	8
Sample Indexes Available:	384
Sequencing Platforms:	Illumina™ NGS

## 1.2 Storage and Stability

- Store the Optics Standard, Optics Solution and pretreatment Solution at -20°C upon arrival to laboratory.
- Store the 384-well Reagent Plate in a sterile and dry environment.

## 1.3 Product Use

- Purified genomic DNA is compatible with this kit.
- Do not use the pretreatment buffer and optics reagents past their expiration date.
- BioQule cartridges cannot be re-used. Take care not to damage or misalign the pipette tips or cartridge tubing. Damaged cartridges or tips may result in assay failure.
- BioQule assays are intended for research use only.
- This manual is a property of Revvity™ Inc.

## 1.4 Warnings and Precautions

We strongly recommend that you read the following warnings and precautions. Periodically, optimizations and revisions are made to the components and manual. Therefore, it is important to follow the most current protocol, which is available on the Revvity website. If you need further assistance, you may contact your local distributor, or contact us at [L3BioQule@Revvity.com](mailto:L3BioQule@Revvity.com).

- Do not use the kit past the expiration date.
- Do not store reagents or pipettes inside the BioQule box.
- Wear gloves and eye protection while setting up the reagent plate for the run.
- Do not place any appendages inside the BioQule while it is running.
- Do not heat Illumina® Unique Dual Index Barcodes above room temperature.
- Do not freeze Illumina® Purification Beads.
- Always vortex the beads to achieve a uniform suspension before pipetting.
- To enable multiplexing, please use the appropriate combination of Illumina® Unique Dual Index Barcodes.
- DNA sample quality may vary between preparations. It is the user's responsibility to utilize high quality DNA. DNA that is heavily nicked or damaged may cause library preparation failure. Absorbance measurements at 260 nm are commonly used to quantify DNA and 260 nm / 280 nm give ratios of 1.8 - 2.0, which usually indicate relatively pure DNA. Other quantification methods using fluorescent dyes may also be used. The user should be aware that contaminating RNA, nucleotides and single-stranded DNA may affect the amount of usable DNA in a sample preparation.

## 1.5 Prior to Starting

- Register your BioQule by sending an email to [L3BioQule@Revvity.com](mailto:L3BioQule@Revvity.com) and get access to training videos, training material, community assay development and software updates.
- Ensure a laboratory temperature of 20° - 25°C (68° - 77°F).
- Identify all reagents and equipment needed during the run in the lab and refrigerators.

## 2 Contents

### 2.1 Kit Contents

The BioQule Illumina® DNA Prep Accessory Kit (P/N 900-000019) contains sufficient materials to prepare 32 libraries. The kit has the following components:

- 4 x BioQule Illumina® DNA Prep Accessory Plates, P/N 810-000020. Each plate comes with a 384 Deep Well Plate, a plate map and a plate loading template insert.
- 4 x BioQule Optics Standard (P/N 820-000060)
- 8 x BioQule Optics Solution (P/N 820-000059)
- 4 x BioQule Pretreatment Buffer (P/N 820-000056)

### 2.2 Additional Equipment, Reagents and Labware

- Equipment
  - BioQule cartridge (Revvity, PN. CLS157064)
  - BioQule NGS Library Prep Instrument (Revvity, PN. CLS155700)
  - Micropipettes: 0.5-10 µl, 2-20 µl, 20-200 µl, 200-1000 µl
  - Microcentrifuge for 0.2 ml tubes
  - Vortexer
  - Plate Centrifuge for SBS Footprint Deep Well Plates
  - Qubit® 2.0, 3.0 or 4.0 Fluoremeter (ThermoFisher Scientific) or other appropriate fluorometer and accessories for quantification of input DNA and final libraries.
  - LabChip GX Touch (Revvity, PN. CLS137031), or equivalent for electrophoretic analysis of nucleic acids.
- Reagents
  - Isopropyl Alcohol (IPA)
  - Nuclease Free Water
  - Illumina® Unique Dual Index Barcodes (Illumina: PN. 20027213, 20027214, 20042666, 20042667)
  - Illumina® DNA Prep Tagmentation Kit (Illumina: PN. 20060059, 96 rxns)
- Supplies and labware
  - Filtered pipette tips, Nuclease Free
  - 0.2 mL PCR strip tubes
  - 10 ml centrifuge tubes

To Order:

- Revvity, [www.revvity.com](http://www.revvity.com)
- Illumina, [www.illumina.com](http://www.illumina.com)
- Fisher Scientific, [www.fishersci.com](http://www.fishersci.com)

## 3 Planning the Run

### 3.1 Workflow and Time Required

BioQule Illumina® DNA Prep Kit for BioQule library prep is a completely automated DNA-Seq library preparation workflow. Each run takes approximately 6 hours with only 30 minutes of hands on time. Figure 1. demonstrates the difference between manual and automated library preparation workflows.

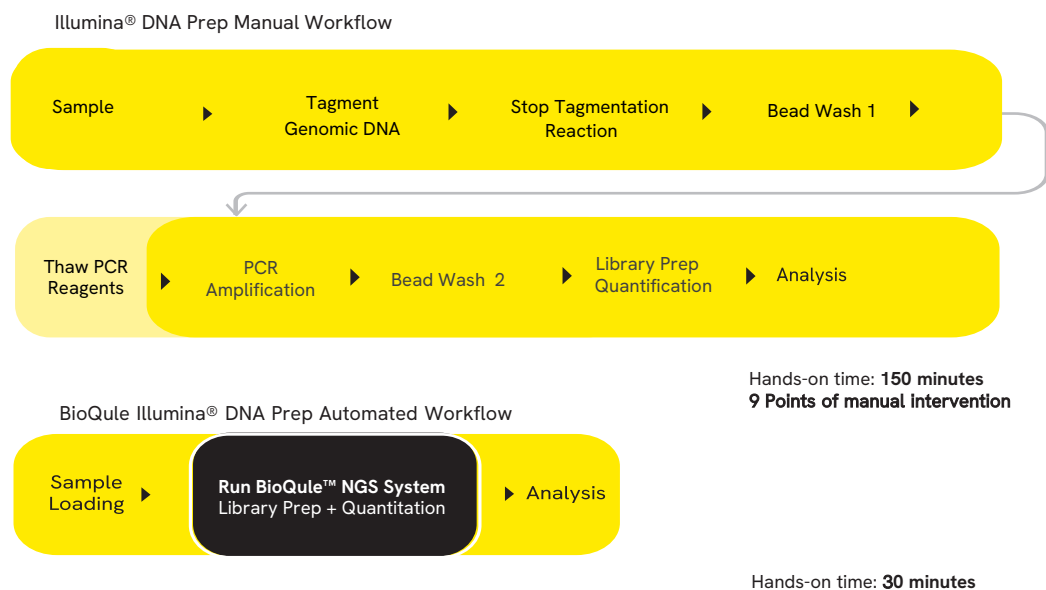


Figure 1. BioQule vs Manual Workflow of Illumina® DNA Prep

### 3.2 Input DNA Requirements

#### DNA Quantity

This kit is compatible with a total DNA input of between 40 ng to 500 ng. Accurate quantification of DNA is required to ensure the minimum input is met. Each set of 8 samples should be normalized to the same input amount to ensure equal amplification for each sample.

#### DNA Purity

DNA Samples must be free of organic solvents (such as phenol, and ethanol), salts used in DNA isolation procedures, any contaminating proteins, and other cellular material. If using a DNA isolation method based on organic solvents, such as TRIzol, we recommend a column purification after isolation. UV absorbance is a common method used for assessing the purity of a DNA sample. The ratio of absorbance at 260 nm to absorbance at 280 nm provides an indication of sample purity. This protocol is optimized for DNA with 260/280 absorbance ratio values of 1.8-2.0, which indicates a pure DNA sample. For a secondary indication of sample purity, use the ratio of absorbance at 260 nm to absorbance at 230 nm. Target a 260/230 ratio of 2.0-2.2

**DNA Integrity**

BioQule Illumina® DNA Prep will generate the best library when DNA samples of high molecular weight and low evidence of degradation are utilized. The BioQule has not been tested with degraded samples. DNA integrity can be determined by utilizing a LabChip GX Touch or equivalent equipment.

**Guidance for Library Amplification**

During the course of run setup, the user will be able to select the number of PCR cycles on the BioQule user interface. This number will default to 8 Cycles but can be adjusted by the user to optimize for specific samples.

### 3.3 Sequencing Recommendations and Guidelines

BioQule Illumina® DNA Prep protocol produces DNA-seq libraries which are compatible with Illumina Sequencing platforms and should be sequenced using the Illumina methodologies following the recommendations for the sequencer being used.

**Index Read Recommendations**

BioQule Illumina® DNA Prep uses 10 base Unique Dual Indexes (UDI) for sample multiplexing. Both Index 1 (i7) and Index 2 (i5) should be sequenced for the detection of "Barcode Hopping". These indexes can be found in Appendix 6.1.

### 3.4 Data Analysis

Once Sequencing data has been generated and parsed, data analysis may be employed according to the requirements of the experiment. If the user requires assistance in this pursuit please contact [L3BioQule@Revvity.com](mailto:L3BioQule@Revvity.com).

### 3.5 Library Storage

Libraries prepared by the BioQule should be stored at a -20°C without a defrost cycle.

## 4 Procedure

### 4.1 Reagent Plate Setup

The steps described below detail how to set up the provided reagent plate for a run on the BioQule. The Plate Loading Guide helps identify which wells will have volume added. The Plate layout seal confirms volumes added into each column.

**Step 1.** Remove the following materials from -20°C storage and thaw for 30 minutes on ice:

- a. Tagmentation Buffer 1 (TB1)
- b. Enhanced PCR Mix (EPM)
- c. Illumina® Index Adaptors
- d. BioQule Optics Standard ( 1 vial)
- e. BioQule Optics Solution ( 2 vials)
- f. BioQule Pretreatment Buffer (1 vial)

**Step 2.** Concurrently, remove the Bead-linked Transposomes (BLT) from 4°C storage and thaw for 30 minutes.

**Step 3.** Prepare the following materials:

- a. New BioQule cartridge
- b. 8 x 15 µl DNA samples in water. Ensure that the DNA concentration is at least 4 ng/µl

*Note: BioQule cartridges cannot be re-used. Take care not to damage or misalign the pipette tips or cartridge tubing. Assay failure may result.*

**Step 4.** Prepare Tagment Stop Buffer (TSB) master mix (8.8x) according to the table below:

Table 1: Tagmentation Stop Buffer Dilution Table

	1x ( µl)	8.8x (µl)
Tagmentation Stop Buffer	13	114.4
Nuclease Free Water	4	35.2
Total	17	149.6

*Note: Vortex to mix. If precipitates are observed, heat at 37°C for 10 minutes, and then vortex until precipitates are dissolved. Mix thoroughly by pipette.*

**Step 5.** Prepare 4mL of 70% IPA. Vortex thoroughly.

Table 2: 70 % IPA

	1x ( µl)	20x (µl)
100 % IPA	140	2800
Nuclease Free Water	60	1200
Total	200	4000



**Step 6.** Prepare Purification Bead Master Mix (8.8x) according to the table below:

Table 3: Purification Bead Mix

	1x ( $\mu$ l)	8.8x ( $\mu$ l)
Illumina® Purification Beads	45	396
Nuclease Free Water	40	352
Total	85	748

*Note: Vortex beads thoroughly before pipetting. Vortex mixture thoroughly until homogenous.*

**Step 7.** Prepare Adapter Mix according to the table below:

Table 4: Adapter Mix

	1x ( $\mu$ l)
Adapter	10
Nuclease free Water	3
Total	13

**Note:**

- Prepare adapter mix separately. DO NOT pool different adapter barcodes.
- Do not reuse barcodes.
- Always take note of the barcodes you have used.
- Immediately put back the adapter plate into  $-20\text{ }^{\circ}\text{C}$  after use.

**Step 8.** For each DNA sample, prepare Tagmentation Master Mix (8.8x) according to the table below:

- Vortex BLT vigorously for 10 seconds to resuspend. Repeat as necessary.
- Vortex TB1 for 10 seconds. Mix thoroughly with a pipette.

Table 5: Tagmentation Master Mix

	1x ( $\mu$ l)
Nuclease Free Water	15
BLT	11
Tagmentation Buffer 1	11
gDNA	15
Total	52

*Note: The 15  $\mu$ l of the gDNA should have a concentration between 2.67 and 33.3 ng/ $\mu$ l which will yield the desired gDNA input amount of 40 - 500 ng. If the gDNA sample is highly concentrated, add a volume lower than 15  $\mu$ l, and add more nuclease free water. If the gDNA concentration is too low, less Nuclease Free Water can be added to allow for adding of a volume greater than 15  $\mu$ l of the gDNA. Total gDNA input amount be between 40 ng and 500 ng.*

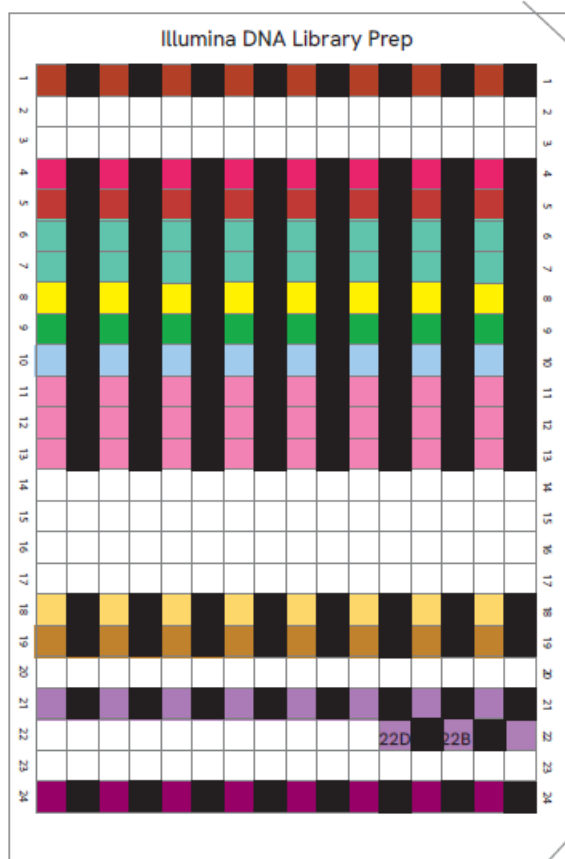


Figure 2. Illumina® DNA Prep Loading Template Guide

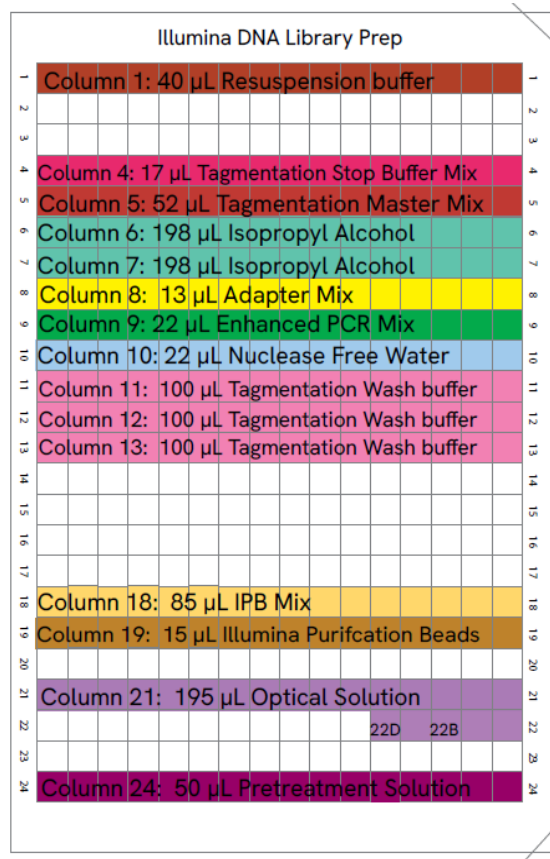


Figure 3. Illumina® DNA Prep Plate Layout

**Step 9.** Follow the plate loading template. Reagents should be loaded in wells marked on the loading template in the following order:

(Do not load the Isopropyl Alcohol until Step 11.)

- Column 1: Load 40 µL of the Resuspension Buffer (RSB).
- Column 4: Load 17 µL Tagment Stop Buffer Mix prepared in Step 4.
- Column 5: Load 52 µL of the Tagmentation Master Mix prepared in Step 8.
- Column 8: Load 13 µL of the Adapter Mix.
- Column 9: Load 22 µL of the EPM.
- Column 10: Load 22 µL of Nuclease free Water (NFW).
- Columns 11, 12 and 13: Vortex and load 100 µL of the Tagmentation Wash Buffer (TWB).
- Column 18: Vortex and load 85 µL of the Purification Bead Mix.
- Column 19: Load 15 µL of undiluted Illumina® purification beads.
- Column 24: Load 50 µL of the Pretreatment Buffer.
- Column 21: Load 195 µL of the Optics Solution.
- Well 22B and 22D: Load 190 µL of the Optics Solution.
- Well 22B: Load 10 µL of the resuspension buffer and pipette mix.
- Well 22D: Load 10 µL of the Optics Standard and pipette mix.

*Note: If large bubbles are present in any column, use a 10 µL pipette tip to pop them gently.*

**Step 10.** Centrifuge the plate for 10 seconds (~1000 rpm)

**Step 11.** After carefully removing the plate from the centrifuge, place it back on the lab bench and load 198  $\mu$ l of 70% IPA into columns 6 and 7. Take extreme care to not spill IPA into adjacent wells, and try not to leave droplets on the plate seal.

*Note: Do not invert or tilt the plate loading template after loading the IPA.*

The Reagent Plate is now ready to run.  
It should look like Figure 4.

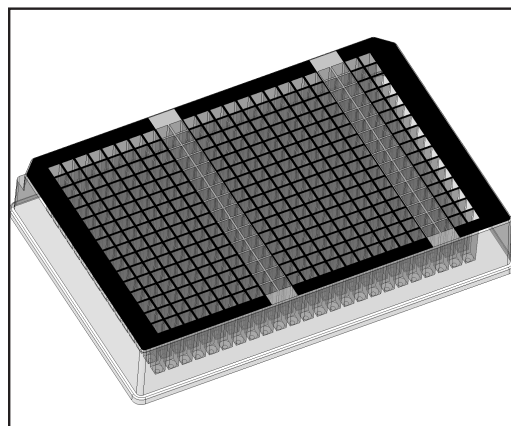


Figure 4. Removing the plate seal from the plate

## 4.2 BioQule Run Setup

The steps described below walk users on how to setup the BioQule instrument and start the run.

**Step 1.** Turn on the BioQule Library Prep System and associated computer, connect the two machines using provided USB cable. Launch the BioQule User Interface on the Computer. Press the Refresh Devices Button to update the Box Connections.

**Step 2.** Select the BioQule machine from the list of options displayed. One computer can run multiple BioQule Boxes. Use the Flash Light button to flash the lights of and identify the selected machine to ensure the correct one is selected. Press Connect to Device to continue.

**Step 3.** On the following screen, select the BioQule Illumina DNA Prep Assay from the list of available assays on the BioQule User Interface.

**Step 4.** Insert the Reagent plate onto the BioQule Plate, as shown in Figure 5.

- a. Make sure the reagent plate is in the correct orientation – the barcode should be facing forwards toward the user, the black seal is up, and the blunt vertex of the 384 well plate should be oriented to the top left.
- b. Press the plate to the left to depress the flat spring on the left side of the x-plate, and then back to depress the flat spring on the back of the x-plate.
- c. Ensure the plate is loaded correctly and is flat against the x-plate

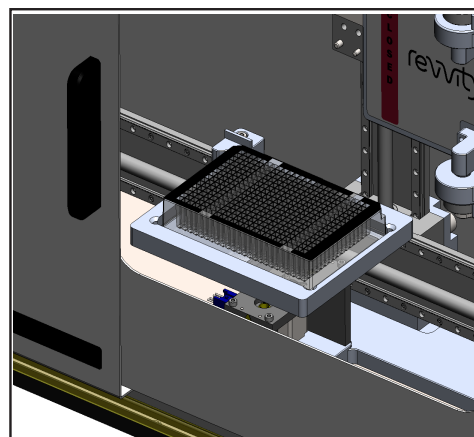


Figure 5. Placing Reagent Plate onto the X-plate

Step 5. The cartridge may now be loaded onto the BioQule instrument.

- a. Begin by removing the PCR door from the instrument.
- b. Hold the cartridge with 2 hands, the cannula array in your left, and the tubing scaffold in your right, make sure the barcode on the tubing scaffold is facing you.
- c. Push the cannula array into the holder, there is an arrow on the pull-tab indication orientation.
- d. Align the eyelets of the cartridge with the 2 posts on the heating element, with one hand on the cartridge at each eyelet, push the cartridge towards the back of the instrument, onto the heating element.

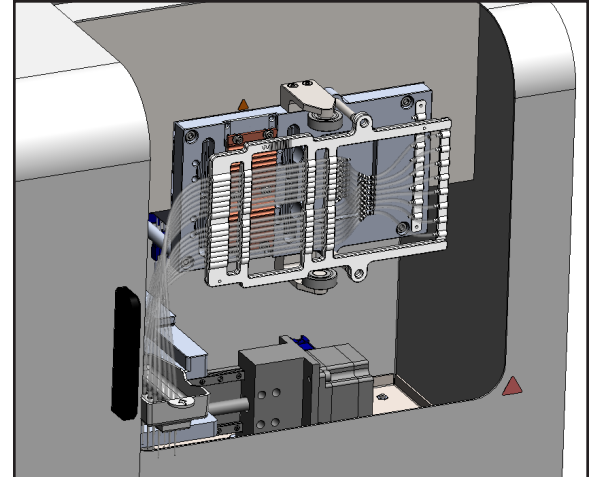


Figure 6. Cartridge insertion into BioQule™

Step 6. Once the cartridge and reagent plate are placed, scan the barcode (using a barcode reader) on each consumable into the correct field on the BioQule UI. Press Enter to confirm Barcodes and then press Next.

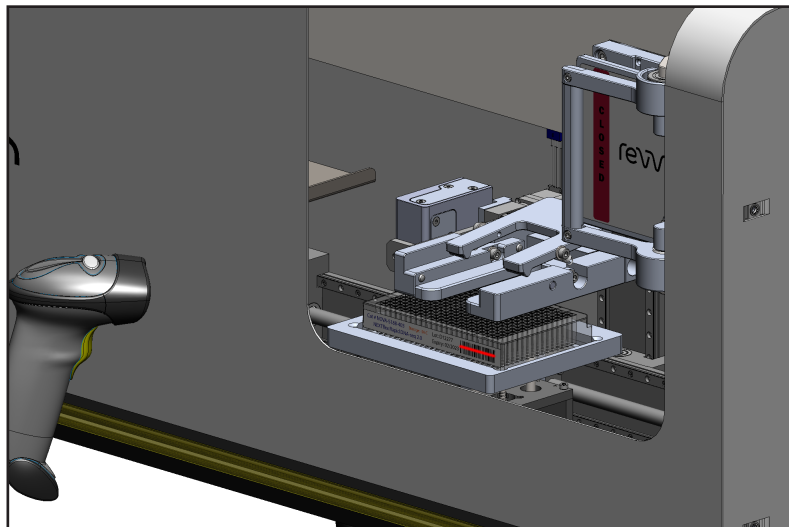


Figure 7. Scan Barcode using Barcode Scanner

Step 7. In the next window, select the ligation time and number of PCR Cycles desired. These values default to 20 minutes and 8 cycles, respectively. The Fragmentation time field is not applicable in this assay.

Step 8. Place the PCR Door with the latch on the left onto BioQule. Then close the PCR door by turning the latch to the right.

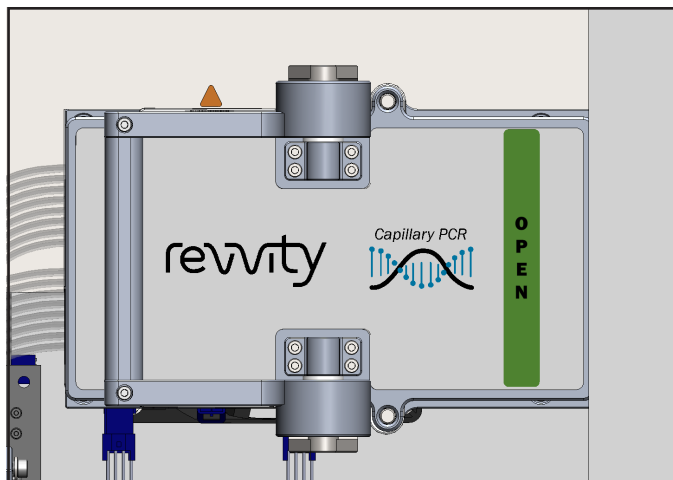


Figure 8. PCR Door Placement

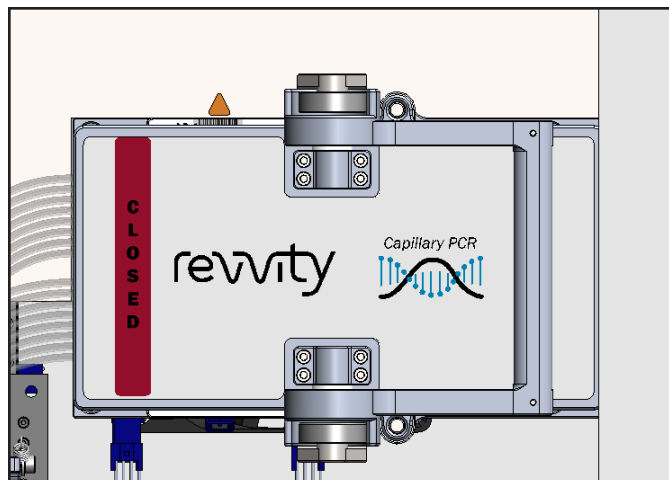


Figure 9. PCR Door Shut

Step 9. Add the names of the Samples and the Sample concentrations to the Spreadsheet. Slide the BioQule door shut and press Run. The Assay will not run unless the door is closed.

- The Assay will now run. It will take approximately 6 hours to complete.
- The Finish button will activate upon completion. DNA Library will be ready and can be found in column 1. The final volume of the library obtained is ~28  $\mu$ L.

Step 11. Spin down the reagent plate and perform quality checks using a LabChip GX Touch prior to sequencing.

*Note: Libraries can be stored in the plate at room temperature for 24 hours. It is recommended to move the libraries to -20°C as soon as reasonably possible but within 24 hours after completion of the protocol.*

Step 12. Assess the quality of the library or pooled libraries using the following method.

- Analyze 1  $\mu$ l library or pooled libraries using the LabChip™\ GX Touch with a High Sensitivity DNA kit.
- The following figures show typical library size profiles with an average fragment size of 550 bp when analyzed.

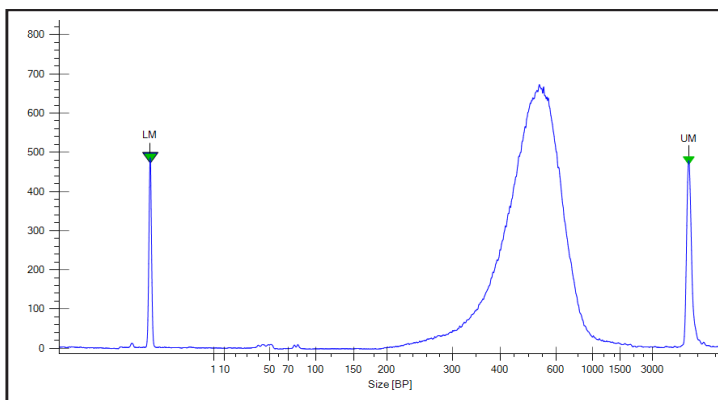


Figure 10. Example Electropherogram Trace for Libraries

*Note: An equivalent electrophoresis platform can be used to analyze the library quality in place of the LabChip GX Touch.*

## 5 Troubleshooting

See Training Videos for explanations of any troubleshooting queries.

### 5.1 Low Volume of Library Generated

Possible causes include:

- Evaporation. Please only leave the plate in the instrument for up to 24 hours. Check the humidity (desire range is 30-50%) and temperature (20 - 25 C) of the lab. The instrument should be far from any devices that vent hot air into the atmosphere.
- Incorrect script was used.
- There may have be bubbles injected into manually loaded wells during the reagent plate loading. Pipetting slowly and into the bottom of the wells is recommended.

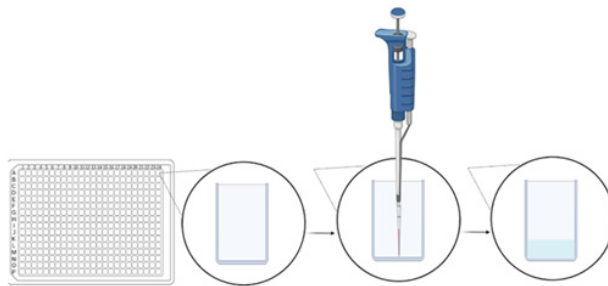


Figure 11. Pipetting directly at bottom of wells

- Large bubbles may have generated in pre-loaded wells upon pipette mixing. Use a 10  $\mu$ l pipette tip to pop them gently.

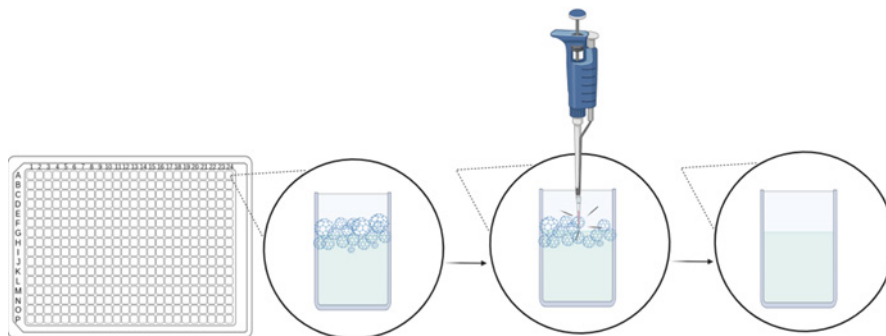


Figure 12. Popping bubbles using a 10  $\mu$ l pipette tip

- BioQule motor stages may have skewed. Please contact [L3BioQule@revvity.com](mailto:L3BioQule@revvity.com).

## 5.2 Low Yield Library Generated

Possible causes include:

- Low quality DNA was used. DNA sample quality may vary between preparations. DNA that is heavily nicked or damaged may cause library preparation failure. Absorbance measurements at 260 nm are commonly used to quantify DNA, and 260 nm / 280 nm ratios of 1.8 - 2.0 usually indicate relatively pure DNA. Other quantification methods using fluorescent dyes may also be used. The user should be aware that contaminating RNA, nucleotides, and single stranded DNA may affect the amount of usable DNA in sample preparation.
- Input DNA amount was incorrect. Input DNA amount should be measured by Qubit or another device. Low input or poor quality of fragmented DNA results in low yield.
- Incorrect pipetting of ligase may result in low yield and high adapters. Ligase is highly viscous. Carefully insert the pipette tip into the bottom of the well, slowly dispense all liquid by pressing the pipette head all the way down, and then move up the pipette tip against the well wall. Once the tip is out of plate well, then depress the pipette.

## 5.3 High Adapter Dimer

Possible causes include:

- Incorrect adapter dilution was used.
- IPA was not freshly made and had expired.
- The Illumina™ Purification Beads were accidentally frozen by mistake or due to a shipment error.
- The Purification Beads were not well resuspended prior to loading onto the BioQule.

# 6 Appendix

## 6.1 Index (UDI) Sequences

Index or Barcode Sequences to be used on Illumina® DNA Prep kits are given below. The actual adapter sequence and what goes onto the sequencing sample sheet is different. The I5 index on the Sample Sheet also differs based on which Sequencer is being used. Use Column 5 if using a NovaSeq 6000 with v1.0 reagent kits, NovaSeq X Series, MiniSeq with Rapid reagents, MiSeq, HiSeq 2000/2500, NextSeq 1000/2000. Use Column 6 if using a iSeq, NovaSeq 6000 with v1.5 reagent kits, MiniSeq, NextSeq 500/550, HiSeq 3000/4000/X, NextSeq 1000/2000.

Table 5: Illumina® UDI Sequences Kit A

Sample Name	I7 Bases in Adapter	I7 Bases on Sample Sheet	I5 Bases in Adapter	I5 Bases on Sample Sheet (1)	I5 Bases on Sample Sheet (2)	Sample Name	I7 Bases in Adapter	I7 Bases on Sample Sheet	I5 Bases in Adapter	I5 Bases on Sample Sheet (1)	I5 Bases on Sample Sheet (2)
UDP0001	CGCTCAGTTC	GAACTGAGCG	TCGTGGAGCG	TCGTGGAGCG	CGTCCACGA	UDP0036	TACGGCCGGT	ACCGGCCGTA	TGGCAATATT	TGGCAATATT	AATATTGCCA
UDP0002	TATCTGACCT	AGGTCAGATA	CTACAAGATA	CTACAAGATA	TATCTTGTAG	UDP0037	GTCCGATTACA	TGTAATCGAC	GATCACCCGCG	GATCACCCGCG	CGCGGTGATC
UDP0003	ATATGAGACG	CGTCTCATAT	TATAGTAGCT	TATAGTAGCT	AGTACTATA	UDP0038	CTGTCTGCAC	GTGCAGACAG	TACCATCCGT	TACCATCCGT	ACGGATGGTA
UDP0004	CTTATGGAAT	ATTCCATAAG	TGCCTGGTGG	TGCCTGGTGG	CCACCAGGCA	UDP0039	CAGCCGATTG	CAATCGGCTG	GCTGTAGGAA	GCTGTAGGAA	TTCTACAGC
UDP0005	TAATCTCGTC	GACGAGATTA	ACATTATCCT	ACATTATCCT	AGGATAATGT	UDP0040	TGACTACATA	TATGTAGTCA	CGCACTAATG	CGCACTAATG	CATTAGTCCG
UDP0006	GCGCGATGTT	AACATCGCGC	GTCCTACTGT	GTCCTACTGT	ACAAGTGGAC	UDP0041	ATTGCCGAGT	ACTCGGCAAT	GACAAGTAA	GACAAGTAA	TTCAAGTTC
UDP0007	AGAGCACTAG	CTAGTGCTCT	TGGAACAGTA	TGGAACAGTA	TACTGTTCCA	UDP0042	GCCATTAGAC	GTCTAATGGC	AGTGGTCAGG	AGTGGTCAGG	CCTGACCACT
UDP0008	TGCCCTTGATC	GATCAAGGCA	CCTTGTTAAT	CCTTGTTAAT	ATTAACAAGG	UDP0043	GGCGAGATGG	CCATCTCGCC	TTCTATGGTT	TTCTATGGTT	AACCATAGAA
UDP0009	CTACTCAGTC	GACTGAGTAG	GTTGATAGTG	GTTGATAGTG	CACTATCAAC	UDP0044	TGGCTCGCAG	CTGCGAGCCA	AATCCGGCCA	AATCCGGCCA	TGGCCGGATT
UDP0010	TCGTCTGACT	AGTCAGACGA	ACCAGCGACA	ACCAGCGACA	TGTCGCTGGT	UDP0045	TAGAATAACG	CGTATTCTA	CCATAAGGTT	CCATAAGGTT	AACCTTATGG
UDP0011	GAACATACGG	CGGTATGTT	CATACACTGT	CATACACTGT	ACAGTGTATG	UDP0046	TAATGGATCT	AGATCCATTA	ATCTCTACCA	ATCTCTACCA	TGGTAGAGAT
UDP0012	CCTATGACTC	GAGTCATAGG	GTGTGGCCT	GTGTGGCCT	AGCGCCACAC	UDP0047	TATCCAGGAC	GTCTGGATA	CGGTGGCGAA	CGGTGGCGAA	TTGCCACCG
UDP0013	TAATGGCAAG	CTTGCCATTA	ATCACAAGG	ATCACAAGG	CCTTCGTGAT	UDP0048	AGTGCCACTG	CAGTGGCACT	TAACAATAGG	TAACAATAGG	CCTATTGTTA
UDP0014	GTGCCGCTTC	GAAGCGCAC	CGGCTCTACT	CGGCTCTACT	AGTAGAGCCG	UDP0049	GTGCAACACT	AGTGTTCAC	CTGGTACACG	CTGGTACACG	CGTGTACCAG
UDP0015	CGGCAATGGA	TCCATTGCCG	GAATGCACGA	GAATGCACGA	TCGTGCATTC	UDP0050	ACATGGTGTG	GACACCATGT	TCAACGTGTA	TCAACGTGTA	TACACGTTGA
UDP0016	GCCGTAACCG	CGGTTACGGC	AAGACTATAG	AAGACTATAG	CTATAGTCTT	UDP0051	GACAGACAGG	CCTGCTGTG	ACTGTTGTA	ACTGTTGTA	TCACAACAGT
UDP0017	AACCATTCTC	GAGAATGGTT	TCGGCAGCAA	TCGGCAGCAA	TTGCTGCCGA	UDP0052	TCTTACATCA	TGATGTAAGA	GTGCGTCTT	GTGCGTCTT	AAGGACGCAC
UDP0018	GGTTCCTCT	AGAGCAACC	CTAATGATGG	CTAATGATGG	CCATCATTAG	UDP0053	TTACAATTC	GGAATTGTA	AGCACATCT	AGCACATCT	AGGATGTCT
UDP0019	CTAATGATGG	CCATCATTAG	GGTTCCTCT	GGTTCCTCT	AGAGGCAACC	UDP0054	AAGCTTATGC	GCATAAGCTT	TTCCGTCGCA	TTCCGTCGCA	TGCGACGGAA
UDP0020	TGGCCCTATC	GATAGGCCGA	CGCACATGGC	CGCACATGGC	GCCATGTGGC	UDP0055	TATTCTCAG	CTGAGGAATA	CTTAACCACT	CTTAACCACT	AGTGGTTAAG
UDP0021	AGTCAACCAT	ATGTTGACT	GGCCTGTCT	GGCCTGTCT	AGGACAGGCC	UDP0056	CTCGTGCTT	AACGCACGAG	GCCTCGGATA	GCCTCGGATA	TATCCGAGGC
UDP0022	GAGCGCAATA	TATTGCCTC	CTGTGTAGG	CTGTGTAGG	CCTAACACAG	UDP0057	TTAGGATAGA	TCTATCTAA	CGTCGACTGG	CGTCGACTGG	CCAGTCGACG
UDP0023	AACAAGCGGT	ACGCCCTGTT	TAAGGAACGT	TAAGGAACGT	ACGTTCTCTA	UDP0058	CCGAAGCGAG	CTCGCTTCGG	TACTAGTCAA	TACTAGTCAA	TTGACTAGTA
UDP0024	GTATGTAGAA	TTCTACATAC	CTAAGTAA	CTAAGTAA	TTACAGTTAG	UDP0059	GGACCAACAG	CTGTTGGTCC	ATAGACGTT	ATAGACGTT	AAGGTTCTAT
UDP0025	TTCTATGGTT	AACCATAGAA	GGCGAGATGG	GGCGAGATGG	CCATCTCGCC	UDP0060	TTCCAGGTAA	TTACCTGGAA	ACAGTCCAG	ACAGTCCAG	CTGGAAGTGT
UDP0026	CCTCGCAACC	GGTTCGAGG	AATAGAGCAA	AATAGAGCAA	TTGCTCTATT	UDP0061	TGATTAGCCA	TGGCTAATCA	AGGCATGTAG	AGGCATGTAG	CTACATGCCT
UDP0027	TGGATGCTTA	TAAGCATCCA	TCAATCCATT	TCAATCCATT	AATGGATTGA	UDP0062	TAACAGTGT	AACACTGTTA	GCAAGTCTCA	GCAAGTCTCA	TGAGACTTGC
UDP0028	ATGCTGTGGT	ACCACGACAT	TCGTATGCGG	TCGTATGCGG	CCGCATACGA	UDP0063	ACCGCGCAAT	ATTGCGCGGT	TTGGCTCCGC	TTGGCTCCGC	GCGGAGCCAA
UDP0029	AGAGTGCAGC	GCCGCACTCT	TCCGACCTCG	TCCGACCTCG	CGAGGTCGGA	UDP0064	GTTCCGCGCA	TGGCGCGAAC	AACTGATACT	AACTGATACT	AGTATCAGTT
UDP0030	TGCCTGGTGG	CCACCAGGCA	CTTATGGAAT	CTTATGGAAT	ATTCCATAAG	UDP0065	AGACACATTA	TAATGTGTCT	GTAAGGCATA	GTAAGGCATA	TATGCCTTAC
UDP0031	TGCGTGTAC	GTGACACGCA	GCTTACGGAC	GCTTACGGAC	GTCGTAAGC	UDP0066	GCGTGGTAT	ATACCAACGC	AATTGTCTGG	AATTGTCTGG	CGCAGCAATT
UDP0032	CATACACTGT	ACAGTGTATG	GAACATACGG	GAACATACGG	CCGTATGTTT	UDP0067	AGCACATCT	AGGATGTGCT	TTACAATTC	TTACAATTC	GGAATTGTAA
UDP0033	CGTATAATCA	TGATTATACG	GTCGATTACA	GTCGATTACA	TGTAATCGAC	UDP0068	TTGTCCGCTG	CACGGAACAA	AACCTAGCAC	AACCTAGCAC	GTGCTAGGTT
UDP0034	TACGCGGCTG	CAGCCGCGTA	ACTAGCCGTG	ACTAGCCGTG	CACGGCTAGT	UDP0069	AAGTACTCCA	TGGAGTACTT	TCTGTGTGGA	TCTGTGTGGA	TCCACACAGA
UDP0035	GCGAGTTACC	GGTAACCTGC	AAGTGTGTA	AAGTGTGTA	TCACCAACTT	UDP0070	ACGTCAATAC	GTATTGACGT	GGAATTCCAA	GGAATTCCAA	TTGGAATTC





Sample Name	17 Bases in Adapter	17 Bases on Sample Sheet	15 Bases in Adapter	15 Bases on Sample Sheet (1)	15 Bases on Sample Sheet (2)
UDP0074	ACCTTATGAA	TTCATAAGGT	TGTTAGAAGG	TGTTAGAAGG	CCTTCTAACA
UDP0075	CGCTGCAGAG	CTCTGCAGCG	GATGGATGTA	GATGGATGTA	TACATCCATC
UDP0076	GTAGAGTCAG	CTGACTCTAC	ACGGCCGTCA	ACGGCCGTCA	TGACGGCCGT
UDP0077	GGATACCAGA	TCTGGTATCC	CGTTGCTTAC	CGTTGCTTAC	GTAAGCAACG
UDP0078	CGCACTAATG	CATTAGTGCG	TGACTACATA	TGACTACATA	TATGTAGTCA
UDP0079	TCCTGACCGT	ACGGTCAGGA	CGGCCTCGTT	CGGCCTCGTT	AACGAGGCCG
UDP0080	CTGGCTTGCC	GGCAAGCCAG	CAAGCATCCG	CAAGCATCCG	CGGATGCTTG
UDP0081	ACCAGCGACA	TGTCGCTGGT	TCGTCTGACT	TCGTCTGACT	AGTCAGACGA
UDP0082	TTGTAACGGT	ACCGTTACAA	CTCATAGCGA	CTCATAGCGA	TCGCTATGAG
UDP0083	GTAAGGCATA	TATGCCTTAC	AGACACATTA	AGACACATTA	TAATGTGTCT
UDP0084	GTCCACTTGT	ACAAGTGGAC	GCGCGATGTT	GCGCGATGTT	AACATCGCGC
UDP0085	TTAGGTACCA	TGGTACCTAA	CATGAGTACT	CATGAGTACT	AGTACTCATG
UDP0086	GGAATTCCAA	TTGGAATTC	ACGTCAATAC	ACGTCAATAC	GTATTGACGT
UDP0087	CATGTAGAGG	CCTCTACATG	GATACCTCCT	GATACCTCCT	AGGAGGTATC
UDP0088	TACACGCTCC	GGAGCGTGTA	ATCCGTAAGT	ATCCGTAAGT	ACTTACGGAT
UDP0089	GCTTACGGAC	GTCCGTAAGC	CGTGTATCTT	CGTGTATCTT	AAGATACACG
UDP0090	CGCTTGAAGT	ACTTCAAGCG	GAACCATGAA	GAACCATGAA	TTCATGGTTC
UDP0091	CGCCTTCTGA	TCAGAAGGCG	GGCCATCATA	GGCCATCATA	TATGATGGCC
UDP0092	ATACCAACGC	GCGTTGGTAT	ACATACTTCC	ACATACTTCC	GGAAGTATGT
UDP0093	CTGGATATGT	ACATATCCAG	TATGTGCAAT	TATGTGCAAT	ATTGCACATA
UDP0094	CAATCTATGA	TCATAGATTG	GATTAAGGTG	GATTAAGGTG	CACCTTAATC
UDP0095	GGTGAATAC	GTATTCCACC	ATGTAGACAA	ATGTAGACAA	TTGTCTACAT
UDP0096	TGGACGGAGG	CCTCCGTCCA	CACATCGGTG	CACATCGGTG	CACCGATGTG

Table 6: Illumina® UDI Sequences Kit B

Sample Name	I7 Bases in Adapter	I7 Bases on Sample Sheet	I5 Bases in Adapter	I5 Bases on Sample Sheet (1)	I5 Bases on Sample Sheet (2)	Sample Name	I7 Bases in Adapter	I7 Bases on Sample Sheet	I5 Bases in Adapter	I5 Bases on Sample Sheet (1)	I5 Bases on Sample Sheet (2)
UDP0097	CTGACCGGCA	TGCCGGTCAG	CCTGATACAA	CCTGATACAA	TTGTATCAGG	UDP0145	GGCCATCATA	TATGATGGCC	GATTGTCATA	GATTGTCATA	TATGACAATC
UDP0098	GAATTGAGTG	CACTCAATTC	TTAAGTTGTG	TTAAGTTGTG	CACAACCTAA	UDP0146	AATTGCTGCG	CGCAGCAATT	ATCCGCTAT	ATCCGCTAT	ATAGCGGAAT
UDP0099	GCGTGTGAGA	TCTCACACGC	CGGACAGTGA	CGGACAGTGA	TCACTGTCCG	UDP0147	TAAGGAACGT	ACGTTCCCTA	GACCGCTGTG	GACCGCTGTG	CACAGCGGTC
UDP0100	TCTCCATTGA	TCAATGGAGA	GCACTACAAC	GCACTACAAC	GTTGTAGTGC	UDP0148	CTATACGGG	CCGCTATAG	TAGGAACCGG	TAGGAACCGG	CCGGTTCCTA
UDP0101	ACATGCATAT	ATATGCATGT	TGGTGCTGG	TGGTGCTGG	CCAGGCACCA	UDP0149	ATTCAGAATC	GATTCTGAAT	AGCGGTGGAC	AGCGGTGGAC	GTCCACCCTG
UDP0102	CAGGCGCCAT	ATGGCGCCTG	TCCACGGCCT	TCCACGGCCT	AGGCGGTGGA	UDP0150	GTAATCTCTA	TAGAGAATAC	TATAGATTCG	TATAGATTCG	CGAATCTATA
UDP0103	ACATAACGGA	TCCGTTATGT	TTGTAGTGT	TTGTAGTGT	TACACTACAA	UDP0151	CCTGATACAA	TTGTATCAGG	ACAGAGGCCA	ACAGAGGCCA	TGGCCTCTGT
UDP0104	TTAATAGACC	GGTCTATTAA	CCACGACACG	CCACGACACG	CGTGTCTGG	UDP0152	GACCGCTGTG	CACAGCGGTC	ATTCCTATTG	ATTCCTATTG	CAATAGGAAT
UDP0105	ACGATTGCTG	CAGCAATCTG	TGTGATGTAT	TGTGATGTAT	ATACATCACA	UDP0153	TTCAGCGTGG	CCACGCTGAA	TATTCTCAG	TATTCTCAG	CTGAGGAATA
UDP0106	TTCTACAGAA	TTCTGTAGAA	GAGCGCAATA	GAGCGCAATA	TATGCGCTC	UDP0154	AACTCCGAAC	GTTCCGAGTT	CGCCTCTGA	CGCCTCTGA	TCAGAAGCGG
UDP0107	TATTGCGTTC	GAACGCAATA	ATCTTACTGT	ATCTTACTGT	ACAGTAAGAT	UDP0155	ATTCGCTAT	ATAGCGGAAT	GCGCAGAGTA	GCGCAGAGTA	TACTCTGCGC
UDP0108	CATGAGTACT	AGTACTCATG	ATGTCGTGGT	ATGTCGTGGT	ACCACGACAT	UDP0156	TGAATATTGC	GCAATATTCA	GGCCCAATT	GGCCCAATT	AATTGGCGCC
UDP0109	TAATTCTACC	GGTAGAATTA	GTAGCCATCA	GTAGCCATCA	TGATGGCTAC	UDP0157	CGCAATCTAG	CTAGATTGCG	AGATATGGCG	AGATATGGCG	CGCCATATCT
UDP0110	ACGCTAATTA	TAATTAGCGT	TGGTTAAGAA	TGGTTAAGAA	TTCTTAACCA	UDP0158	AACCGCATCG	CGATGCGGTT	CCTGCTTGGT	CCTGCTTGGT	ACCAAGCAGG
UDP0111	CCTGTTAAT	ATTAACAAGG	TGTTGTTCTG	TGTTGTTCTG	ACGAACAACA	UDP0159	CTAGTCCGGA	TCCGACTAG	GACGAACAAT	GACGAACAAT	ATTGTTCTGC
UDP0112	GTAGCCATCA	TGATGGCTAC	CCAACAACAT	CCAACAACAT	ATGTTGTTGG	UDP0160	GCTCCGTCAC	GTGACGGAGC	TGGCGGTCCA	TGGCGGTCCA	TGGACCGCCA
UDP0113	CTTGTAATTC	GAATTACAAG	ACCGGCTCAG	ACCGGCTCAG	CTGAGCCGGT	UDP0161	AGATGGAATT	AATCCATCT	CTTCAGTTAC	CTTCAGTTAC	GTAAGTGAAG
UDP0114	TCCAATCTA	TAGAATTGGA	GTTAATCTGA	GTTAATCTGA	TCAGATTAAC	UDP0162	ACACCGTAA	TTAACGGTGT	TCCTGACCGT	TCCTGACCGT	ACGGTCAGGA
UDP0115	AGAGCTGCCT	AGGCAGCTCT	CGGCTAACGT	CGGCTAACGT	ACGTTAGCCG	UDP0163	GATAACAAGT	ACTTGTATC	CGCGCTAGA	CGCGCTAGA	TGTAGCGCGG
UDP0116	CTTCGCCGAT	ATCGCGGAAG	TCCAAGAATT	TCCAAGAATT	AATTCTTGA	UDP0164	CTGTACACG	CGGTACCAG	AGGATAAGTT	AGGATAAGTT	AACTTATCT
UDP0117	TCGGTCACGG	CCGTGACCGA	CCGAACGTTG	CCGAACGTTG	CAACGTTCCG	UDP0165	CGAAGGTTAA	TTAACCTTCG	AGGCCAGACA	AGGCCAGACA	TGCTGGCCT
UDP0118	GAACAAGTAT	ATACTGTTC	TAACCGCGA	TAACCGCGA	TCGCGGGTTA	UDP0166	ATCGCATATG	CATATGCAT	CCTTGAACGG	CCTTGAACGG	CCGTTCAAGG
UDP0119	AATTGGCGGA	TCCGCCAATT	CTCCGTGCTG	CTCCGTGCTG	CAGCAGGAG	UDP0167	ATCATAGGCT	AGCCTATGAT	CACCACCTAC	CACCACCTAC	GTAGGTGGTG
UDP0120	GGCCTGTCT	AGGACAGGCC	CATTCCAGCT	CATTCCAGCT	AGCTGGAATG	UDP0168	GATTGTCATA	TATGACAATC	TTGCTTGTAT	TTGCTTGTAT	ATACAAGCAA
UDP0121	TAGGTTCTCT	AGAGAACCTA	GGTTATGCTA	GGTTATGCTA	TAGCATAACC	UDP0169	CCAACAACAT	ATGTTGTTGG	CAATCTATGA	CAATCTATGA	TCATAGATTG
UDP0122	ACACAATATC	GATATTGTGT	ACCACACGGT	ACCACACGGT	ACCGTGTGGT	UDP0170	TTGTTGGTGC	GCACCACCAA	TGGTACTGAT	TGGTACTGAT	ATCAGTACCA
UDP0123	TTCTGTGACG	CGTACAGGAA	TAGGTTCTCT	TAGGTTCTCT	AGAGAACCTA	UDP0171	GCGAAGCGCT	AGCGTTCGCG	TTATCCAAC	TTATCCAAC	GTTGGATGAA
UDP0124	GGTAACGACG	CTGCGTTACC	TATGGCTCGA	TATGGCTCGA	TCGAGCCATA	UDP0172	CAACCGGAGG	CCTCCGTTG	CATAACCCA	CATAACCCA	TGGTGTATG
UDP0125	TCCACGGCCT	AGGCGTGGGA	CTCGTGCGTT	CTCGTGCGTT	AACGACGAG	UDP0173	AGCGGTGGAC	GTCCACCCTG	TCCTATTAGC	TCCTATTAGC	GCTAATAGGA
UDP0126	GATACCTCCT	AGGAGGTATC	CCAGTTGGCA	CCAGTTGGCA	TGCCAACTGG	UDP0174	GACGAACAAT	ATTGTTGCTC	TCTCTAGATT	TCTCTAGATT	AATCTAGAGA
UDP0127	CAACGTCAGC	GCTGACGTTG	TGTTGCAATT	TGTTGCAATT	AATGCGAACA	UDP0175	CCACTGGTCC	GGACCAGTGG	CGCGAGCCTA	CGCGAGCCTA	TAGGCTCGCG
UDP0128	CGGTTATTAG	CTAATAACCG	AACCGCATCG	AACCGCATCG	CGATGCGGTT	UDP0176	TGTTAGAAGG	CCTTCTAACA	GATAAGCTCT	GATAAGCTCT	AGAGCTTATC
UDP0129	CGCGCTAGA	TCTAGGCGCG	CGAAGGTTAA	CGAAGGTTAA	TTAACCCTCG	UDP0177	TATATTGAG	CTCGAATATA	GAGATGTCGA	GAGATGTCGA	TCGACATCTC
UDP0130	CTTTGGCTAT	ATAGCCAAGA	AGTGCCACTG	AGTGCCACTG	CAGTGGCACT	UDP0178	CGCGACGATC	GATGTCGCG	CTGGATATGT	CTGGATATGT	ACATATCCAG
UDP0131	TCACACGAA	TTGCGTGTGA	GAACAAGTAT	GAACAAGTAT	ATACTGTTC	UDP0179	GCCTCGATA	TATCCGAGGC	GGCCAATAAG	GGCCAATAAG	CTTATTGGCC
UDP0132	AACGTTACAT	ATGTAACGTT	ACGATTGCTG	ACGATTGCTG	CAGCAATCGT	UDP0180	TGAGACAGCG	CGCTGTCTCA	ATTACTCACC	ATTACTCACC	GGTGAATAT
UDP0133	CGGCCTCGTT	AACGAGGCCG	ATACCTGGAT	ATACCTGGAT	ATCCAGGTAT	UDP0181	TGTTGCGATT	AATGCGAACA	AATTGGCGGA	AATTGGCGGA	TCCGCCAATT
UDP0134	CATAACACCA	TGTTGTTATG	TCCAATTCTA	TCCAATTCTA	TAGAATTGGA	UDP0182	TCCAAGAATT	AATTCTTGA	TTGTCAACT	TTGTCAACT	AAGTTGACAA
UDP0135	ACAGAGGCCA	TGCGCTCTGT	TGAGACAGCG	TGAGACAGCG	CGCTGTCTCA	UDP0183	GCTGTAGGAA	TTCTACAGC	GGCGAATTCT	GGCGAATTCT	AGAATTCGCC
UDP0136	TGGTGCTGG	CCAGGCACCA	ACGCTAATTA	ACGCTAATTA	TAATTAGCGT	UDP0184	ATACCTGGAT	ATCCAGGTAT	CAACGTCAGC	CAACGTCAGC	GCTGACGTTG
UDP0137	TAGGAACCGG	CCGGTTCCTA	TATATTGAG	TATATTGAG	CTCGAATATA	UDP0185	GTTGGACCGT	ACGGTCCAAC	TCTTACATCA	TCTTACATCA	TGATGTAAGA
UDP0138	AATATTGGCC	GGCCAATATT	CGGTCGATA	CGGTCGATA	TATCGGACCG	UDP0186	ACCAAGTTAC	GTAATTGGT	CGCCATACCT	CGCCATACCT	AGGTATGGCG
UDP0139	ATAGGTATTC	GAATACCTAT	ACAATAGAGT	ACAATAGAGT	ACTCTATTGT	UDP0187	GTGTGGCGCT	AGCGCCACAC	CTAATGTCTT	CTAATGTCTT	AAGACATTAG
UDP0140	CCTTCACGTA	TACGTGAAGG	CGGTTATTAG	CGGTTATTAG	CTAATAACCG	UDP0188	GGCAGTAGCA	TGCTACTGCC	CAACCGGAGG	CAACCGGAGG	CCTCCGGTTG
UDP0141	GGCCAATAAG	CTTATTGGCC	GATAACAAGT	GATAACAAGT	ACTTGTATT	UDP0189	TGCGGTGTTG	CAACACCGCA	GGCAGTAGCA	GGCAGTAGCA	TGCTACTGCC
UDP0142	CAGTAGTTGT	ACAACACTGT	AGTTATCACA	AGTTATCACA	TGTGATAACT	UDP0190	GATTAAGGTG	CACCTTAATC	TAGGATAGCA	TAGGATAGCA	CTATCTCTAA
UDP0143	TTCATCCAAC	GTTGATGAA	TTCCAGGTAA	TTCCAGGTAA	TTACCTGGAA	UDP0191	CAACATTCAA	TTGAATGTTG	CGCAATCTAG	CGCAATCTAG	CTAGATTGCG
UDP0144	CAATTGGATT	AATCCAATTG	CATGTAGAGG	CATGTAGAGG	CCTCTACATG	UDP0192	GTGTTACCGG	CCGGTAACAC	GAGTTGACT	GAGTTGACT	AGTACAACCT



Table 7: Illumina® UDI Sequences Kit C

Sample Name	17 Bases in Adapter	17 Bases on Sample Sheet	15 Bases in Adapter	15 Bases on Sample Sheet (1)	15 Bases on Sample Sheet (2)
UDP0193	TATCATGAGA	TCTCATGATA	AACACGTGGA	AACACGTGGA	TCCACGTGTT
UDP0194	CTTGGCCTCG	CGAGGCCAAG	GTGTTACCGG	GTGTTACCGG	CCGGAACAC
UDP0195	GTCTCTGAA	TTACAGAGAC	AGATTGTTAC	AGATTGTTAC	GTAACAATCT
UDP0196	CCATCCACGC	CGGTGGATGG	TTGACCAATG	TTGACCAATG	CATTGGTCAA
UDP0197	ACAACCAGGA	TCCTGGTGTG	CTGACCGGCA	CTGACCGGCA	TGCCGGTCAG
UDP0198	AGCAGAATTA	TAATCTGCT	TCTCATCAAT	TCTCATCAAT	ATTGATGAGA
UDP0199	CAGTCGTGCG	CGCACGACTG	GGACCAACAG	GGACCAACAG	CTGTGGTGTC
UDP0200	GTCTAACCTC	GAGGTAGAC	AATGTATTGC	AATGTATTGC	GCAATACATT
UDP0201	GAACTCGGTT	AACCGAGTTC	GATCTCTGGA	GATCTCTGGA	TCCAGAGATC
UDP0202	AGTTATCACA	TGTGATAACT	CAGGCGCCAT	CAGGCGCCAT	ATGGCGCCTG
UDP0203	GTAGCATACT	AGTATGCTAC	TTAATAGACC	TTAATAGACC	GGTCTATTAA
UDP0204	CTTCAGTTAC	GTAACCTGAAG	GGAGTCGCGA	GGAGTCGCGA	TCGCGACTCC
UDP0205	AGTCCGAGGA	TCCTGGGACT	AACGCCAGAG	AACGCCAGAG	CTCTGGCGTT
UDP0206	ACAGTTCCAG	CTGGAACCTG	CGTAATTAAC	CGTAATTAAC	GTTAATTACG
UDP0207	CCGCATATTC	GAATATGCGG	ACGAGACTGA	ACGAGACTGA	TCAGTCTCGT
UDP0208	TTATCCGATC	GATCGGATAA	GTATCGGCCG	GTATCGGCCG	CGGCCGATAC
UDP0209	ATAGTCTAGC	GCTAGACTAT	AATACGACAT	AATACGACAT	ATGTCGTATT
UDP0210	TATAGTAGCT	AGCTACTATA	GTTATATGGC	GTTATATGGC	GCCATATAAC
UDP0211	ACTCCGGTGG	CCACCGGAGT	GCCTGCCATG	GCCTGCCATG	CATGGCAGGC
UDP0212	GTGCGTAAG	CTTACCACAC	TAAGACCTAT	TAAGACCTAT	ATAGGCTTAA
UDP0213	GATATCCTAA	TTAGGATATC	TATACCATGG	TATACCATGG	CCATGGTATA
UDP0214	TCGCGTATAA	TTATACGCGA	GCCGTCTGTT	GCCGTCTGTT	AACAGACGGC
UDP0215	ATTCTAAGCG	CGCTTAGAAT	CAGAGTGATA	CAGAGTGATA	TATCACTCTG
UDP0216	AGCGCTTCGG	CCGAAGCGCT	TGCTAACTAT	TGCTAACTAT	ATAGTTAGCA
UDP0217	GTTGATAGTG	CACTATCAAC	TCAGTTAATG	TCAGTTAATG	CATTAACTGA
UDP0218	AATAGAGCAA	TTGCTCTATT	GTGACCTTGA	GTGACCTTGA	TCAAGGTCAC
UDP0219	CTAACTGTAA	TTACAGTTAG	ACATGCATAT	ACATGCATAT	ATATGCATGT
UDP0220	GCGTACTTAG	CTAAGTACGC	AACATACCTA	AACATACCTA	TAGGTATGTT
UDP0221	TACCGAACTA	TAGTTCGGTA	CCATGTGTAG	CCATGTGTAG	CTACACATGG
UDP0222	GTAGTAATAG	CTATTACTAC	GAGTCTCTCC	GAGTCTCTCC	GGAGAGACTC
UDP0223	GGTTATGCTA	TAGCATAACC	GCTATGCGCA	GCTATGCGCA	TGCGCATAGC
UDP0224	ACAATAGAGT	ACTCTATTGT	ATCGCATATG	ATCGCATATG	CATATGCGAT
UDP0225	GCCTCCACTA	TAGTGAAGCG	AGTACCTATA	AGTACCTATA	TATAGTACT
UDP0226	AGATATGGCG	CGCCATATCT	GACCGGAGAT	GACCGGAGAT	ATCTCCGGTC
UDP0227	AATATGAAGC	GCTTCATATT	CGTTCAGCCT	CGTTCAGCCT	AGGCTGAACG
UDP0228	TAGCGCTAGT	ACTAGCGCTA	TTACTTCTCT	TTACTTCTCT	GAGGAAGTAA
UDP0229	AGTTAAGAGC	GCTCTTAACT	CACGTCCACC	CACGTCCACC	GGTGAGCAGT
UDP0230	CAGATACCAC	GTGGTATCTG	GCTACTATCT	GCTACTATCT	AGATAGTAGC
UDP0231	ACGGCCGTCA	TGACGGCCGT	AGTCAACCAT	AGTCAACCAT	ATGGTTGACT
UDP0232	GTAATTAAGT	CAGTAATTAAC	CGAGGCGGTA	CGAGGCGGTA	TACCGCCTCG
UDP0233	AAGTCTTGTA	TACAAGACTT	CAGGTGTTCA	CAGGTGTTCA	TGAACACCTG
UDP0234	GTCACCACAG	CTGTGGTGAC	GACAGACAGG	GACAGACAGG	CCTGTCTGTC
UDP0235	ATTAGTGGAG	CTCCACTAAT	TGTAATTGTT	TGTAATTGTT	AACAAGTACA
UDP0236	TGCTAACTAT	ATAGTTAGCA	CTCTAAGTAG	CTCTAAGTAG	CTACTTAGAG
UDP0237	TAAGACCTAT	ATAGTCTTAA	GTCACCACAG	GTCACCACAG	CTGTGGTGAC
UDP0238	TGGTTAAGAA	TTCTTAACCA	TCTACATACC	TCTACATACC	GGTATGAGA
UDP0239	ACTTCTCTT	AAGGAAGAGT	CACGTTAGGC	CACGTTAGGC	GCCTAACGTG
UDP0240	GTCTCTTCC	GGAAGGAGAC	TGGTGAGTCT	TGGTGAGTCT	AGACTCACCA
UDP0241	TCCCGTTC	TGAACGCGGA	CTTCGAAGGA	CTTCGAAGGA	TCCTCGAAG

Sample Name	17 Bases in Adapter	17 Bases on Sample Sheet	15 Bases in Adapter	15 Bases on Sample Sheet (1)	15 Bases on Sample Sheet (2)
UDP0242	AGTTGCAGG	CCTGCAACCT	GTAGAGTCAG	GTAGAGTCAG	CTGACTCTAC
UDP0243	GAACCATGAA	TTATCGTTC	GACATTGTCA	GACATTGTCA	TGACAATGTC
UDP0244	TTGAGAGGAT	ATCCTCTCAA	TCCGCAAGGC	TCCGCAAGGC	GCCTTCCGGA
UDP0245	TGGTCTAGTG	CACTAGACCA	ACTGCCTTAT	ACTGCCTTAT	ATAAGGCAGT
UDP0246	AGTGGATAAT	ATTATCCACT	TACGCACGTA	TACGCACGTA	TACGTGCGTA
UDP0247	GGCACGCCAT	ATGGCGTGCC	CGCTTGAAGT	CGCTTGAAGT	ACTTCAAGCG
UDP0248	GATCTCTGGA	TCCAGAGATC	CTGCACCTCA	CTGCACCTCA	TGAAGTGCAG
UDP0249	TGCTGGACAT	ATGTCACGCA	CAGCGGACAA	CAGCGGACAA	TTGTCCGCTG
UDP0250	CCGAACGTTG	CAACGTTCCG	GGATCCGCAT	GGATCCGCAT	ATGCGGATCC
UDP0251	ATTAATACGC	CGTATTAAT	TGCGGTGTTG	TGCGGTGTTG	CAACACCGCA
UDP0252V2	CCAGATTCGG	CCGAATCTGG	ATGAATCAAG	ATGAATCAAG	CTTGATTCT
UDP0252	TAGTCAACA	GTTGTGACTA	ACATAACGGA	ACATAACGGA	TCCGTTATGT
UDP0253	GGTATTGAGA	TCTCAATACC	GACGTTCCGC	GACGTTCCGC	CGCGAACGTC
UDP0254	CAAGATGCTT	AAGCATCTTG	CATTCAACAA	CATTCAACAA	TTGTGAATG
UDP0255	ACGAGACTGA	TCAGTCTCGT	CACGGATTAT	CACGGATTAT	ATAATCCGTC
UDP0256	TTATCTTGCA	TGCAAGATAA	TTGAGGACGG	TTGAGGACGG	CCGCTCTCAA
UDP0257	AGATTGTTAC	GTAACAATCT	CTCTGTATAC	CTCTGTATAC	GTATACAGAG
UDP0258V2	TATACCATGG	CCATGGTATA	TCTCGCGGAG	TCTCGCGGAG	CTCCGCGAGA
UDP0258	TGTACCCTG	CAGCGGTAGA	GCAACAGGTG	GCAACAGGTG	CACCTGTTGC
UDP0259	AACGGTATGA	TCATACCGTT	GGTAACGCG	GGTAACGCG	CTGCGTTACC
UDP0260	CAATGGCGCC	GGCGCCATTG	ACCGCGCAAT	ACCGCGCAAT	ATTGCGCGGT
UDP0261	CTAATTCGCT	AGCGAATAG	AGCCGGAACA	AGCCGGAACA	TGTTCCGCT
UDP0262	CATGGTCTAA	TTAGACCATG	TCCTAGGAAG	TCCTAGGAAG	CTTCTAGGA
UDP0263	ATACTGTGTG	CACACAGTAT	TTGAGCCTAA	TTGAGCCTAA	TTAGGCTCAA
UDP0264	GCCGACAAGA	TCTTGTCCGC	CCACCTGTGT	CCACCTGTGT	ACACAGGTGG
UDP0265	CGAGGCGGTA	TACCGCCTCG	CCTCGCAACC	CCTCGCAACC	GGTTGCGAGG
UDP0266	GATATAACAG	CTGTTATATC	GTATAGCTGT	GTATAGCTGT	ACAGCTATAC
UDP0267	TCGCGGTTA	TAACCGGCGA	GCTACATTAG	GCTACATTAG	CTAATGATAG
UDP0268	AGACTCTCTT	AAGAGAGTCT	TACGAATCTT	TACGAATCTT	AAGATTCGTA
UDP0269	GCTCGCTTAC	GTAGGCGGAG	TAGGAGCGCA	TAGGAGCGCA	TGCGCTCTTA
UDP0270	AGGATAAGTT	AACTTATCTT	GTACTGGCGT	GTACTGGCGT	ACGCGAGTAC
UDP0271	GAGACATAAT	ATTATGTCTC	AGTTAAGAGC	AGTTAAGAGC	GCTCTTAACT
UDP0272	AGCTGTATATA	TATAACAGCT	TCGCGTATAA	TCGCGTATAA	TTATACGCGA
UDP0273	GTATCATGG	CCAATGATAC	GAGTGTGCG	GAGTGTGCG	CGGCACACTC
UDP0274	AATAGCCTC	GAGGCTTATT	CTAGTCCGGA	CTAGTCCGGA	TCCGGACTAG
UDP0275	CCGCTTAGCT	AGCTAAGCGG	ATTAATACGC	ATTAATACGC	GCGTATTAAT
UDP0276	TCCTAGGAAG	CTTCTAGGTA	CCTAGAGTAT	CCTAGAGTAT	ATACTTAGG
UDP0277	TCACAGATCG	CGATCTGTGA	TAGGAAGACT	TAGGAAGACT	AGTCTTCTTA
UDP0278	ACTTGCCAC	GTGACAAGT	CCGTGGCCTT	CCGTGGCCTT	AAGGCCACGG
UDP0279	TGTACTTGT	AACAAGTACA	GGATATATCC	GGATATATCC	GGATATATCC
UDP0280	CACCTAATCT	AGATAAGTG	CACCTCTTGG	CACCTCTTGG	CCAAGAGGTG
UDP0281	CAGAGTGATA	TATCACTCTG	AACGTTACAT	AACGTTACAT	ATGTAACGTT
UDP0282	GGCGAATTCT	AGAATTCGCC	CGGCAAGCTC	CGGCAAGCTC	GAGCTTGCCG
UDP0283	AGTGGTCAGG	CCTGACCACT	TCTTGGCTAT	TCTTGGCTAT	ATAGCCAAGA
UDP0284	CATTCCAGCT	AGCTGGAATG	ACGGAATGCG	ACGGAATGCG	CGCATCCGT
UDP0285	CTCGTTATCA	TGATAACGAG	GTCCGCGAGG	GTCCGCGAGG	CCTGCGGAAC
UDP0286	CCTTACTATG	CATAGTAAGG	ACCAAGTTAC	ACCAAGTTAC	GTAACCTGGT
UDP0287	AGAAGCCAAT	ATTGGCTTCT	TGGCTCGCAG	TGGCTCGCAG	CTGCGAGCCA
UDP0288	TAATCGGATC	GTACCGATTA	AACTAACGTT	AACTAACGTT	AACGTTAGTT



Table 8: Illumina® UDI Sequences Kit D

Sample Name	17 Bases in Adapter	17 Bases on Sample Sheet	15 Bases in Adapter	15 Bases on Sample Sheet (1)	15 Bases on Sample Sheet (2)	Sample Name	17 Bases in Adapter	17 Bases on Sample Sheet	15 Bases in Adapter	15 Bases on Sample Sheet (1)	15 Bases on Sample Sheet (2)
UDP0289V2	GCTACTATCT	AGATAGTAGC	GGCACGCCAT	GGCACGCCAT	ATGGCGTGCC	UDP0335	CCGGAATCAT	ATGATTCCGG	CTCTGACGTG	CTCTGACGTG	CACGTCAGAG
UDP0289	GGAATTGTTT	GAACAATTCC	TAGAGTTGGA	TAGAGTTGGA	TCCAACCTTA	UDP0336	TTGAGCCTAA	TTAGGCTCAA	TCGAATGGAA	TCGAATGGAA	TTCCATTGGA
UDP0290V2	GTCTTCTAAT	ATTAGAAGAC	GCAGGCTGGA	GCAGGCTGGA	TCCAGCCTGC	UDP0337	CCACCTTACA	TGTAAGGTGG	AAGGCCTTGG	AAGGCCTTGG	CCAAGGCCTT
UDP0290	CCGGACCACA	TGTGGTCCGG	AGAGCACTAG	AGAGCACTAG	CTAGTGCTCT	UDP0338	GTTCAGTTG	CAACTGCAAC	TGAACGCAAC	TGAACGCAAC	GTTCGCTTCA
UDP0291V2	ATGTGCGAGC	GCTCGCATAT	ATGGCTTAAT	ATGGCTTAAT	ATTAAGCCAT	UDP0339	TCACCTAGT	ACATGAGTGA	CCGCTTAGCT	CCGCTTAGCT	AGCTAAGCGG
UDP0291	GACTTAGAAG	CTTCTAAGTC	ACTCTACAGG	ACTCTACAGG	CCTGTAGAGT	UDP0340	GACTGGTTGC	GCAACCAGTC	CACCGAGGAA	CACCGAGGAA	TTCTCGGTG
UDP0292	TGGCAATATT	AATATTGCCA	CGGTGACACC	CGGTGACACC	GGTGTACCCG	UDP0341	ATCGTCGCTC	GAGCGACGAT	CGTATAATCA	CGTATAATCA	TGATTATACG
UDP0293	GAATGCACGA	TCGTGCATTC	CGCTTGGTAT	CGCTTGGTAT	ATACCAACGC	UDP0342	GGTGCCTTCG	CGAACGCACC	ATGACAGAAC	ATGACAGAAC	GTTCGTGCAT
UDP0294	CGTGTATCTT	AAGATACACG	TGTGCTAACA	TGTGCTAACA	TGTTAGCACACA	UDP0343	CGGCGTAAGA	TCTTACGCCG	ATTCATTGCA	ATTCATTGCA	TGCAATGAAT
UDP0295	ATTCATTGCA	TGCAATGAAT	CCAGAAGTAA	CCAGAAGTAA	TTACTTCTGG	UDP0344	GACATCAGCT	AGCTGATGTC	TCATGCTCTG	TCATGCTCTG	CAGGACATGA
UDP0296	TCCTTCATAG	CTATGAAGGA	CTTATACCTG	CTTATACCTG	CAGTATAAG	UDP0345	ACTAATTCAG	CTGAATTAGT	AATTCGATCG	AATTCGATCG	CGATCGAATT
UDP0297	TCTAGTCTTC	GAAGACTAGA	ACTAGAACTT	ACTAGAACTT	AAGTCTAGT	UDP0346	TTCTCCTCTA	TAAGGAGGAA	TTCCGACATT	TTCCGACATT	AATGTCGGAA
UDP0298	CTCGACTCCT	AGGAGTCGAG	TTAGGCTTAC	TTAGGCTTAC	GTAAGCCTAA	UDP0347	TGTGTAAGCT	AGCTTACACA	TGGCAGCACC	TGGCAGCACC	GGTGTGCCA
UDP0299	AGTGAGTGAA	TTCACTCACT	TATCATGAGA	TATCATGAGA	TCTCATGATA	UDP0348	GTGGCTGTTT	AACCAGCCAC	GCCACAGCAC	GCCACAGCAC	GTGCTGTGGC
UDP0300	GAAGCGGACC	GGTCCGCTTC	CTCACACAAG	CTCACACAAG	CTTGTGTGAG	UDP0349	TCGACTTAAG	CCTAAGTCGA	CAGTAGTTGT	CAGTAGTTGT	ACAACACTG
UDP0301V2	CAAGCCACTA	TAGTGGCTTG	AGTTACTTGG	AGTTACTTGG	CCAAGTAACT	UDP0350	CACGTTAGGC	GCCTAACGTG	AGCTCTCAAG	AGCTCTCAAG	CTTGAGAGCT
UDP0301	GCTCTCGTTG	CAACGAGAGC	GAATTGAGTG	GAATTGAGTG	CACTCAATTC	UDP0351	TGAAGTAAGT	ACTACTTCA	TCTGGAATTA	TCTGGAATTA	TAATCCAGA
UDP0302	GGACCTCAAT	ATTGAGGTCC	CGGATTATAT	CGGATTATAT	ATATAATCCG	UDP0352	ACGGAATGCG	CGCATTCCGT	ATTAGTGGAG	ATTAGTGGAG	CTCCACTAAT
UDP0303	GAGTCTCTCC	GGAGAGACTC	TTGAAGCAGA	TTGAAGCAGA	TCTGCTTCAA	UDP0353	GTGTGATATC	GATATCACAC	GACTATATGT	GACTATATGT	ACATATAGTC
UDP0304	AACGAGCGG	CCGCTCCGTT	TACGCGAAG	TACGCGAAG	CTTCGCCGTA	UDP0354	ACACAGCGCT	AGCGTGTGT	CGTTCGGAAC	CGTTCGGAAC	GTTCGGAACG
UDP0305	TGTGATGAT	ATACATCACACA	TCTCCATTGA	TCTCCATTGA	TCAATGGAGA	UDP0355	AGCGCGGTGA	TCACCGCGCT	TCGATACTAG	TCGATACTAG	CTAGTATCGA
UDP0306	AACATACCTA	TAGGTATGTT	CGAGACCAAG	CGAGACCAAG	CTTGGTCTCG	UDP0356	CAAGGCTATC	GATAGCCTTG	TACCACAATG	TACCACAATG	CATTGTGGTA
UDP0307	GTGCTAGGTG	CACCTAGCAC	TGCTGGACAT	TGCTGGACAT	ATGTCACGA	UDP0357	TGCCTCCAGG	CCTGGACGCA	TGGTATACCA	TGGTATACCA	TGGTATACCA
UDP0308	CATACTTGAA	TTCAAGTATG	GATGGTATCG	GATGGTATCG	CGATACCATC	UDP0358	AGGTGCGTAA	TTACGCACCT	GCTCTCGTTG	GCTCTCGTTG	CAACGAGAGC
UDP0309	CTTGTCTTAA	TTAAGACAAG	GGCTTAATTG	GGCTTAATTG	CAATTAAGCC	UDP0359	GCAGCAACGA	TCGTTGCTGC	GTCTCGTGAA	GTCTCGTGAA	TTCACGAGAC
UDP0310	AAGAGAGGTG	CACCTCTCTT	CTCGACTCCT	CTCGACTCCT	AGGAGTCGAG	UDP0360	ATCCTTGTGC	CGACAAGGAT	AAGGCCACCT	AAGGCCACCT	AGGTGGCCTT
UDP0311	TGCACAGAGAA	TTCTCGTGCA	ATACACAGAG	ATACACAGAG	CTCTGTGTAT	UDP0361	GAAGGTACAC	GTGTACCTTC	CTGTGAGCTA	CTGTGAGCTA	TAGCTCACAG
UDP0312	ACTTCTAGC	GCTAGGAAGT	TCTCGGACGA	TCTCGGACGA	TCGTCCGAGA	UDP0362	TTGGCCAGGT	ACCTGGCCAA	TCACAGATCG	TCACAGATCG	CGATCTGTGA
UDP0313	GTGCTATTA	TTAATAGCAC	ACCACGCTCG	ACCACGCTCG	CAGACGTGGT	UDP0363	AGGCAGACA	TGCTGCGCT	AGAAGCCAAT	AGAAGCCAAT	ATTGGCTTCT
UDP0314	AGCGTGAATG	CATTACGCT	GTTGACTCA	GTTGACTCA	TGAGTACAAC	UDP0364	AGCATTAACT	AGTTAATGCT	ACTGCAGCCG	ACTGCAGCCG	CGGCTGCAGT
UDP0315	CCTTAGTGCC	GGCACTAAGG	TCAGGTCAAC	TCAGGTCAAC	GTTGACCTGA	UDP0365	ATTACTCACC	GGTGAGTAAT	AACATCTAGT	AACATCTAGT	ACTAGATGTT
UDP0316	TGTACCGAAT	ATTCGGTACA	AGTCCGAGGA	AGTCCGAGGA	TCCTCGGACT	UDP0366	GCGCAGAGTA	TACTCTGCGC	CCTTACTATG	CCTTACTATG	CATAGTAAGG
UDP0317	GGAGATTAGT	ACTAATCTCC	CACCTAATCT	CACCTAATCT	AGATTAAGTG	UDP0367	CGCCATACCT	AGGTATGGCG	GTGGCGAGAC	GTGGCGAGAC	GTCTCGCCAC
UDP0318	TACTAACACA	TGTGTTAGTA	TACTCTGTTA	TACTCTGTTA	TAACAGAGTA	UDP0368	GCAGCTGGA	TCCAGCTGCG	GCCAGATCCA	GCCAGATCCA	TGGATCTGGC
UDP0319	TAGTCTGTTG	CAACGACCTA	GCGACTCGAT	GCGACTCGAT	ATCGAGTCGC	UDP0369	GTTATATGCG	GCCATATAAC	ACACAATATC	ACACAATATC	GATATTGTGT
UDP0320	ATGCCAGCCG	CGGTCCGAT	CTAGGCAAGG	CTAGGCAAGG	CCTTGCCCTAG	UDP0370	CACCTCGACT	AGTGCAGGTG	TGGAGGTAAT	TGGAGGTAAT	ATTACCTCCA
UDP0321	CTAGCGTCGA	TCGACGCTAG	CCTCTTGGAA	CCTCTTGGAA	TTCGAAGAGG	UDP0371	ACCGGCTCAG	CTGAGCCGGT	CCTTACAGTA	CCTTACAGTA	TACGTGAAGG
UDP0322	TGCCTACGAG	CTCGTAGGCA	TCATCCTCTT	TCATCCTCTT	AAGAGGATGA	UDP0372	ATAGACCGTT	AACGGTCTAT	CTATACCGGG	CTATACCGGG	CCGCGTATAG
UDP0323	ACTAGAAGCTT	AAGTTCTAGT	GGTAAGATAA	GGTAAGATAA	TTATCTTACC	UDP0373	TGAACGCAAC	GTTCGCTTCA	GTTCGAGTTG	GTTCGAGTTG	CAACTGCAAC
UDP0324	CACCTCTTGG	CCAAGAGGTG	AACGAGCCAG	AACGAGCCAG	CTGGCTCGTT	UDP0374	GTGGTTGAAG	CTTCAACCAC	TTATGCGCTT	TTATGCGCTT	AGGCGCATAA
UDP0325	AAGCAGATAT	ATATCTGCTT	TAGACAATCT	TAGACAATCT	AGATTGTCTA	UDP0375	ACTGAATAGA	TCTATTGAGT	TCTCAGTACA	TCTCAGTACA	TGACTGAGAG
UDP0326	GCCAGATCCA	TGGATCTGGC	CAATGCTGAA	CAATGCTGAA	TTCAGCATTG	UDP0376	GGACGCTCTG	CAAGACGTCC	AGTATACGGA	AGTATACGGA	TCCGTATACT
UDP0327	TTGATTCAA	TTGAATCCAA	GTCACGGTGT	GTCACGGTGT	ACACCGTGAC	UDP0377	GTTGACTCA	TGAGTACAAC	ACGCTTGGAC	ACGCTTGGAC	GTCCAAGCGT
UDP0328	ACTAGCCGTG	CACGGCTAGT	GGGTACAAG	GGGTACAAG	CTGTACACC	UDP0378	AGAACCOCGG	CCGCGTCTCT	GGAGTAGATT	GGAGTAGATT	AATCTACTCC
UDP0329	CGGCAAGCTC	GAGCTTCCCG	AGGTTGCAGG	AGGTTGCAGG	CCTGCAACCT	UDP0379	CAGTATCAAT	ATTGATACTG	TACACGCTCC	TACACGCTCC	GGAGCGTGTGA
UDP0330	GAAGCTAGCT	AGCTAGCTTC	TAATACGGAG	TAATACGGAG	CTCCGTATTA	UDP0380	TCCATAATCC	GGATTATGGA	TCCGATAGAG	TCCGATAGAG	CTCTATCGGA
UDP0331	ACAAGGATTG	CAATCCTTGT	CGAAGACGCA	CGAAGACGCA	TGCGTCTTCG	UDP0381	ATGAGAACCA	TGGTCTCAT	CTCAAGGCCG	CTCAAGGCCG	CGGCCTGAG
UDP0332	GCAACAGGTG	CACCTGTTGC	ATTGACACAT	ATTGACACAT	ATGTGTCAAT	UDP0382	TCGTGTTTGA	TCAACCACGA	CAAGTTCATA	CAAGTTCATA	TATGAACCTG
UDP0333	CAAGGTGACG	CGTCACTTGG	CAGCCGATTG	CAGCCGATTG	CAATCGGCTG	UDP0383	CAAGTTCATA	TATGAACCTG	AATCCTTAGG	AATCCTTAGG	CCTAAGGATT
UDP0334	ACCAGTCATT	AATGACTGGT	TCTCACGCGT	TCTCACGCGT	ACGCGTGAGA	UDP0384	CTTAACCACT	AGTGGTTAAG	GGTGAATAC	GGTGAATAC	GTATTCCACC



## 7 Technical Assistance

For help with any of our products, please contact Revvity Technical Support at +1 203-925-4602 (direct) or 800.762.4000 (toll-free, U.S. only) or email [L3BioQule@revvity.com](mailto:L3BioQule@revvity.com), or fill out a Customer Support form on our website [www.revvity.com/customer-support](http://www.revvity.com/customer-support).

## 8 Revision History

Date	Revision	Notes
July 19th 2024	E	Updated Electropherograms

The Revvity logo is displayed in a lowercase, sans-serif font.