

Takara Bio USA

# ThruPLEX® DNA-Seq HV User Manual

Cat. Nos. R400740, R400741  
(031320)

---

**Takara Bio USA, Inc.**

1290 Terra Bella Avenue, Mountain View, CA 94043, USA

U.S. Technical Support: [techUS@takarabio.com](mailto:techUS@takarabio.com)

United States/Canada  
800.662.2566

Asia Pacific  
+1.650.919.7300

Europe  
+33.(0)1.3904.6880

Japan  
+81.(0)77.565.6999

## Table of Contents

|  |    |
|--|----|
| I. Introduction.....   | 3  |
| A. Overview.....   | 3  |
| B. Principle.....  | 4  |
| C. ThruPLEX DNA-Seq HV Workflow.....                               | 5  |
| II. List of Components.....  | 6  |
| A. Components.....   | 6  |
| B. Shipping and Storage Conditions.....                            | 6  |
| C. Additional Materials Required.....                              | 6  |
| D. Optional Materials.....   | 7  |
| III. General Considerations.....                                   | 7  |
| A. Sample Requirements.....  | 7  |
| B. General Recommendations.....                                    | 9  |
| C. Safety Guidelines.....  | 11 |
| IV. Protocols.....   | 12 |
| A. Protocol: ThruPLEX DNA-Seq HV Library Preparation.....          | 12 |
| B. Library Processing for Illumina Next-Generation Sequencing..... | 16 |
| Appendix A. Indexing Reagents.....                                 | 19 |
| A. Overview.....   | 19 |
| B. Components.....   | 19 |
| C. ThruPLEX HV Unique Dual Index Sequences.....                    | 19 |
| Appendix B. Troubleshooting Guide.....                             | 22 |

## Table of Figures

|   |    |
|---|----|
| Figure 1. ThruPLEX DNA-Seq HV single-tube library preparation workflow.....   | 3  |
| Figure 2. ThruPLEX DNA-Seq HV technology uses a three-step, single-tube reaction that starts with fragmented double-stranded DNA or cfDNA (5 ng to 200 ng)..... | 4  |
| Figure 3. Overview of ThruPLEX DNA-Seq HV library preparation for Illumina NGS, starting with fragmented DNA....  | 5  |
| Figure 4. Fragment size distribution of cfDNA isolated from plasma.....   | 9  |
| Figure 5. Example of real-time analysis of library amplification using ThruPLEX DNA-Seq HV.....   | 10 |
| Figure 6. Bioanalyzer analysis of libraries prepared using ThruPLEX DNA-Seq HV.....   | 18 |

## Table of Tables

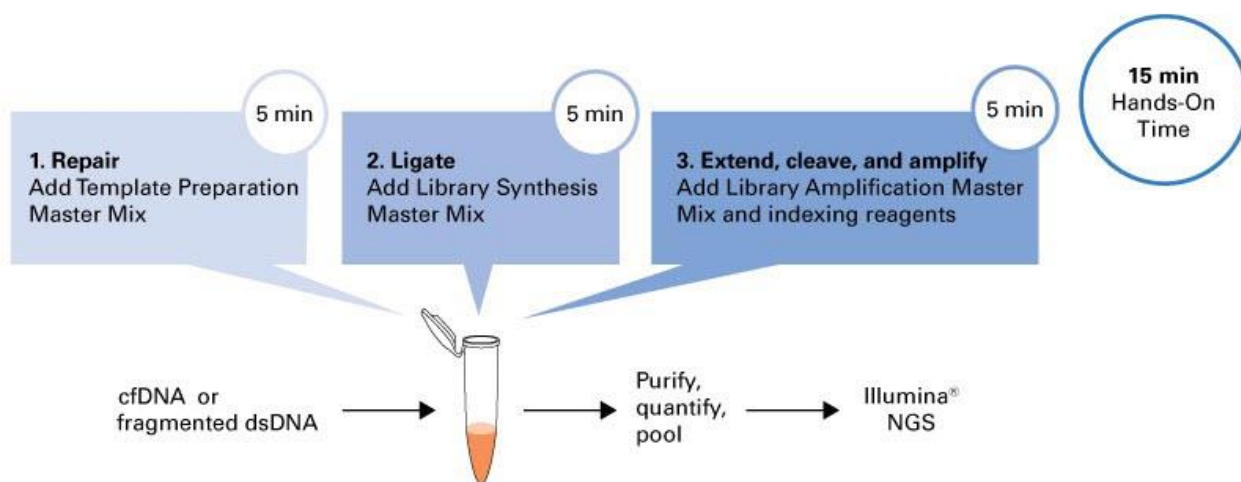
|   |    |
|---|----|
| Table I. ThruPLEX DNA-Seq HV Kit Contents.....                                  | 6  |
| Table II. DNA Sample Requirements.....  | 7  |
| Table III. Recommended DNA Purification Kits.....                               | 7  |
| Table IV. Estimated gene copies based on input amount and allele frequency..... | 8  |
| Table V. ThruPLEX HV UDI - Set A plate layout.....                              | 19 |
| Table VI. ThruPLEX HV UDI sequences.....  | 20 |
| Table VII. Troubleshooting Guide for the ThruPLEX Tag-Seq HV Kits.....          | 22 |

## I. Introduction

### A. Overview

The ThruPLEX HV DNA-Seq HV kit is designed to provide up to 96 indexed libraries for higher multiplexing capabilities on Illumina® NGS platforms. ThruPLEX DNA-Seq HV chemistry is engineered and optimized to generate DNA libraries with high molecular complexity and balanced GC representation from input volumes of up to 30 µl. Inputs of 5 ng to 200 ng of fragmented double-stranded DNA are required for library preparation. The entire three-step workflow takes place in a single tube or well, in about two hours (Figure 1). No intermediate purification steps and no sample transfers are necessary, thus preventing handling errors and loss of valuable samples. With high library diversity, ThruPLEX DNA-Seq HV libraries excel when combined with target enrichment and deliver high-quality sequencing results.

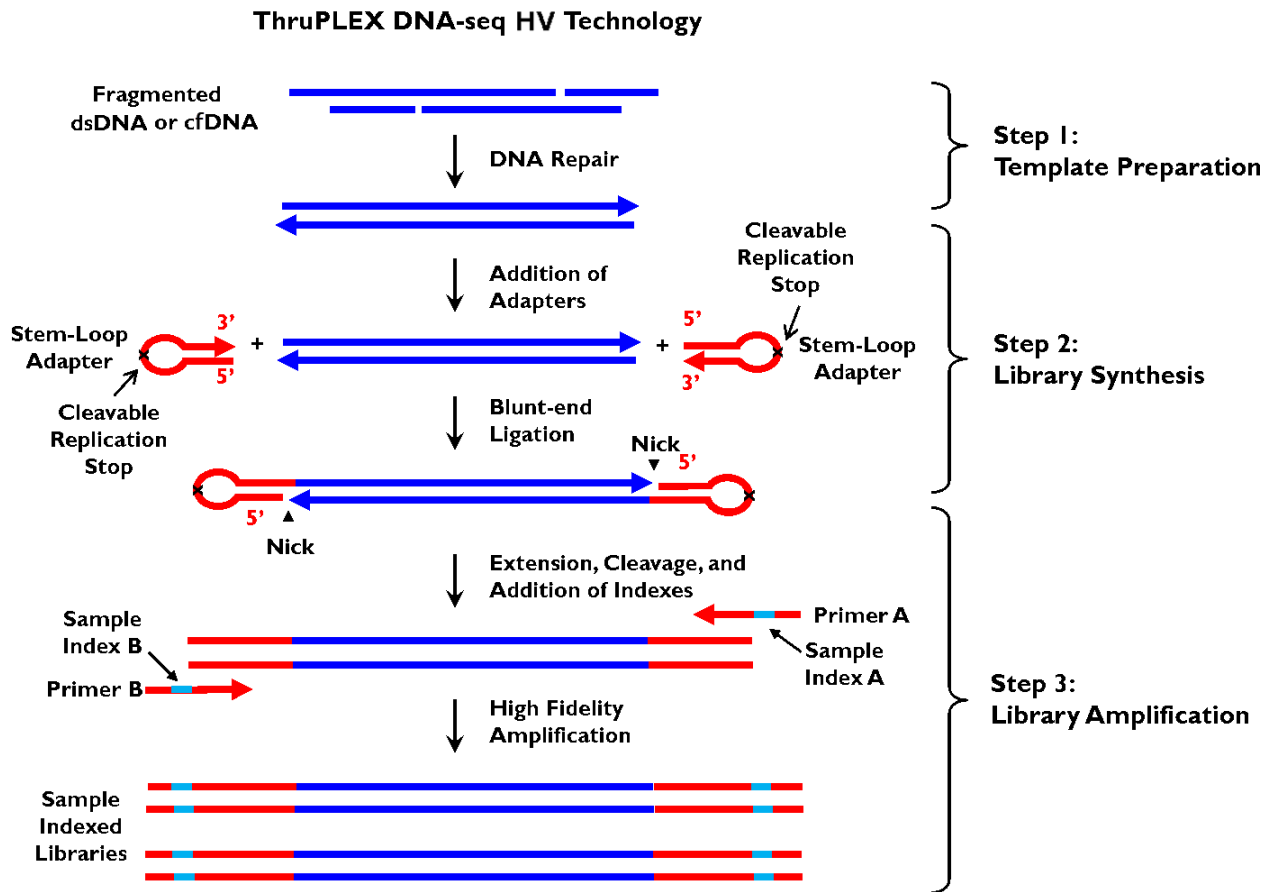
Pairing ThruPLEX DNA-Seq HV with ThruPLEX HV unique dual indexes (UDIs) adds the capability of multiplexing up to 96 NGS-ready libraries. Once purified and quantified, the resulting libraries are ready for Illumina NGS instruments using standard Illumina sequencing reagents and protocols. The kit provides excellent results for high-coverage deep sequencing, such as *de novo* sequencing, whole genome resequencing, whole exome sequencing, and/or other enrichment techniques. It is ideally suited for use with small fragments of DNA such as cell-free plasma DNA or damaged DNA from formalin-fixed paraffin-embedded (FFPE) tissue.



**Figure 1. ThruPLEX DNA-Seq HV single-tube library preparation workflow.** The ThruPLEX DNA-Seq HV workflow consists of three simple steps that take place in the same PCR tube or well and eliminates the need to purify and transfer the sample material.

**B. Principle**

The ThruPLEX DNA-Seq HV kit is based on our patented ThruPLEX HV technology (Figure 2). Unlike other NGS library preparation kits, which are based on ligation of Y-adapters, ThruPLEX HV uses stem-loop adapters to construct high-quality libraries in a fast and efficient workflow. In the first step, Template Preparation, the DNA is repaired and yields molecules with blunt ends. In the next step, Library Synthesis, stem-loop adapters with blocked 5' ends are ligated with high efficiency to the 5' end of the genomic DNA, leaving a nick at the 3' end. The adapters cannot ligate to each other and do not have single-strand tails, both of which contribute to nonspecific background found with many other NGS preparations. In the final step, the 3' ends of the genomic DNA are extended to complete library synthesis, and Illumina-compatible indexes are added through high-fidelity amplification. Any remaining free adapters are destroyed. Hands-on time and risk of contamination are minimized by using a single tube and eliminating intermediate purifications.



**Figure 2.** ThruPLEX DNA-Seq HV technology uses a three-step, single-tube reaction that starts with fragmented double-stranded DNA or cfDNA (5 ng to 200 ng). Stem-loop adapters are blunt-end ligated to repaired input DNA. These molecules are extended, then amplified to include barcodes using a high-fidelity polymerase to yield an indexed Illumina NGS library.

## C. ThruPLEX DNA-Seq HV Workflow

The ThruPLEX DNA-Seq HV workflow is highly streamlined (Figure 3) and consists of the following three steps:

- **Template Preparation** for efficient repair of the fragmented double-stranded DNA input.
- **Library Synthesis** for ligation of our patented stem-loop adapters.
- **Library Amplification** for extension of the template, cleavage of the stem-loop adapters, and amplification of the library. Illumina-compatible indexes are also introduced using a high-fidelity, highly-processive, low-bias DNA polymerase.

The three-step ThruPLEX DNA-Seq HV workflow takes place in a single tube or well and is completed in about two hours.

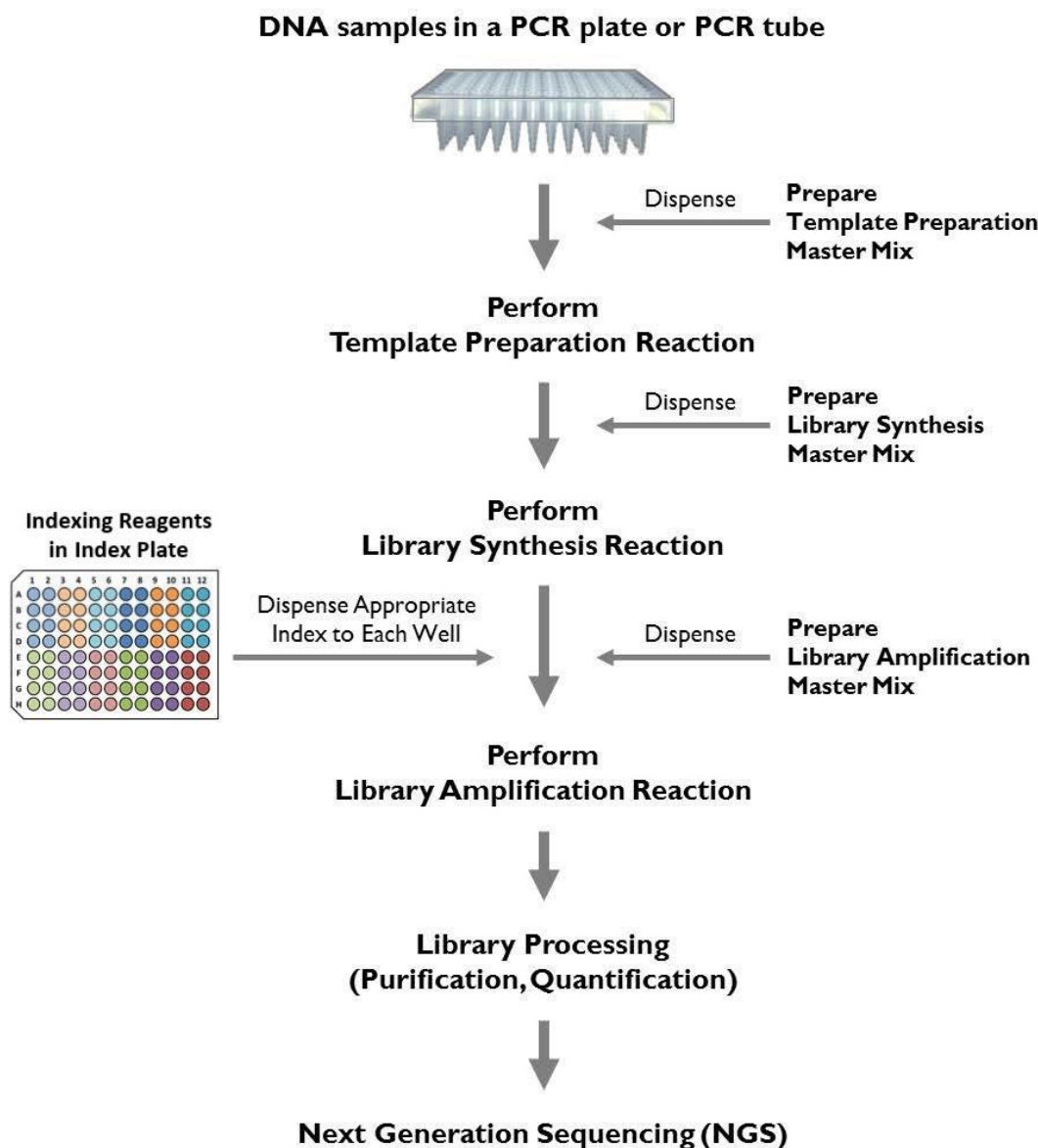


Figure 3. Overview of ThruPLEX DNA-Seq HV library preparation for Illumina NGS, starting with fragmented DNA.

## II. List of Components

### A. Components

Table I. ThruPLEX DNA-Seq HV Kit Contents

| Name                                    | Cap color  | Storage | R400741<br>(24 rxns)         | R400740<br>(96 rxns)         |
|---|------------|---------|------------------------------|------------------------------|
| PBD1                                    | Blue       | -20°C   | 1 tube                       | 1 tube                       |
| PED1                                    | Blue       | -20°C   | 1 tube                       | 1 tube                       |
| SBD1                                    | White      | -20°C   | 1 tube                       | 1 tube                       |
| SED1                                    | White      | -20°C   | 1 tube                       | 1 tube                       |
| ABD1                                    | Amber tube | -20°C   | 1 tube                       | 4 tubes                      |
| AED1                                    | Violet     | -20°C   | 1 tube                       | 1 tube                       |
| Control Fragmented Human gDNA (5 ng/μl) | N/A        | -20°C   | 1 tube                       | 1 tube                       |
| Nuclease-Free Water                     | Clear      | -20°C   | 1 tube                       | 1 tube                       |
| ThruPLEX HV UDI*                        |            | -20°C   | 1 Dual Index<br>Plate (24 D) | 1 Dual Index<br>Plate (96 D) |

\*included in bundle part numbers R400740 & R400741; also sold separately as R400738 & R400739

### B. Shipping and Storage Conditions

ThruPLEX DNA-Seq HV is shipped on dry ice. The kit should be stored at -20°C upon arrival.

### C. Additional Materials Required

- ThruPLEX HV Indexing module (part number R400738 or R400739) if ThruPLEX DNA-Seq HV was purchased as a core components kit
- Hot-lid PCR thermal cycler (real-time instrument optional)

**NOTE:** See Thermal Cycler Considerations in Section III.B.1.

- Centrifuge
  - PCR tubes or 96-well nuclease-free thin-wall PCR plates
- NOTE:** Select appropriate tubes or plates that are compatible with the thermal cyclers and/or real-time thermal cyclers used. Use appropriate caps or sealing films and seal thoroughly to eliminate evaporation during cycling conditions. Evaporation could reduce robustness and reproducibility of the reactions.
- 1.5 ml low adhesion microcentrifuge tubes
  - PCR plate seals (if using plates)
  - Single-channel pipette: 10 μl, 20 μl, and 200 μl
  - Multi-channel pipettes: 20 μl and 200 μl
  - Low-binding filter pipette tips: 10 μl, 20 μl, 200 μl
  - Low-binding aerosol barrier tips
  - Low TE (10 mM Tris, 0.1 mM EDTA, pH 8.0)
  - 80% (v/v) ethanol: freshly made for each experiment
  - Magnetic Separator, such as SMARTer-Seq™ Magnetic Separator - PCR Strip (Takara Bio, Cat. No. 635011)
  - Fluorometer, such as Qubit, for library quantification
  - Agencourt AMPure XP beads (Beckman Coulter, Cat. No. A63880)

## ThruPLEX® DNA-Seq HV User Manual

**NOTE:** Agencourt AMPure XP beads need to come to room temperature before the container is opened. Therefore, we strongly recommend aliquoting the beads upon receipt, and then refrigerating the aliquots. Individual tubes can be removed for each experiment, allowing them to come to room temperature more quickly (~30 min). This aliquoting process is also essential for minimizing the chances of bead contamination.

Immediately prior to use, vortex the beads until they are well dispersed. The color of the liquid should appear homogeneous. Confirm that there is no remaining pellet of beads at the bottom of the tube. Mix well to disperse before adding the beads to your reactions. The beads are viscous, so pipette them slowly.

### D. Optional Materials

- 20X fluorescent dye, such as EvaGreen, in water (Biotium, Cat. No. 31000-T)
- Calibration Dye (such as Bio-Rad Laboratories, Cat. No. 170-8780)
- Reference Dye (if required by real time instrument)
- qPCR-based library quantification kit for Illumina NGS libraries: Library Quantification Kit (Takara Bio, Cat. No. 638324)
- Bioanalyzer or TapeStation, for library size distribution

## III. General Considerations

### A. Sample Requirements

Table II. DNA Sample Requirements

| DNA Sample Requirements |  |
|-------------------------|--|
| Nucleic Acid Source     | Fragmented double-stranded DNA or cfDNA<br>Cells, plasma, urine, other biofluids, FFPE, tissues, fresh tissues, frozen tissues |
| Type                    | Mechanically sheared; enzymatically fragmented; low-molecular-weight cell-free DNA   |
| Molecular Size          | <1,000 bp  |
| Input Amount            | 5 ng to 200 ng   |
| Input Volume            | 30 $\mu$ l   |
| Input Buffer            | $\leq$ 10 mM Tris, $\leq$ 0.1 mM EDTA  |

#### 1. General Guidelines

DNA samples must be fragmented dsDNA in order to be used with ThruPLEX DNA-Seq HV. Fragmented double-stranded DNA (gDNA), degraded DNA from sources such as FFPE, cfDNA from plasma, or other biofluids are suitable. This kit is not for use with single-stranded DNA (ssDNA) or RNA.

#### 2. DNA Isolation

The table below lists recommended kits for isolation of common sample types. For additional recommendations, please contact Takara Bio Technical Support.

Table III. Recommended DNA Purification Kits.

| Sample type                                | Recommended kit           | Catalog Nos.                        |
|--|---------------------------|-------------------------------------|
| FFPE tissue                                | NucleoSpin DNA<br>FFPE XS | 740980.10, 740980.50,<br>740980.250 |
| Plasma, urine, etc.                        | NucleoSnap DNA<br>Plasma  | 740300.10, 740300.50                |
| Mammalian cells and tissues                | NucleoSpin Tissue         | 740952.10, 740952.50,<br>740952.250 |
| Mammalian cells and tissues<br>(low input) | NucleoSpin Tissue<br>XS   | 740901.10, 740901.50,<br>740901.250 |

**3. Input DNA Amount**

The recommended input amount is 5 ng to 200 ng of DNA quantified by Qubit Fluorometer or equivalent methods. When working with cfDNA, quantification of the mononucleosomal cfDNA fragments by Bioanalyzer run is recommended. Use an appropriate input amount of DNA to ensure sufficient variant copies are available for the library preparation process to achieve the desirable detection sensitivity. In general, detection of alleles present at low frequencies requires higher input amount of DNA.

**Table IV. Estimated gene copies based on input amount and allele frequency.**

| Estimated genome copies available for library preparation |                              |  |     |      |
|---|------------------------------|--|-----|------|
| Input amount  | Total haploid genome copies* | Total variant copies at indicated allele frequency |     |      |
|   |                              | 5%   | 1%  | 0.5% |
| 100 ng  | 33,333                       | 1,666  | 333 | 166  |
| 50 ng   | 16,666                       | 833  | 166 | 83   |
| 10 ng   | 3,333                        | 166  | 33  | 16   |
| 5 ng  | 1,666                        | 83   | 16  | 8    |

\*Calculated using 3 pg as the mass of a haploid genome. The genomic complexity of plasma samples is highly variable. All numbers are rounded down to the nearest whole number.

**4. Input Volume**

The maximum input sample volume is 30 µl. If a sample is in a larger volume, the DNA must be concentrated into 30 µl or less. Care should be taken to ensure the buffer concentration is appropriate (see below).

**5. Input Buffer**

Input DNA must be eluted or re-suspended in a low-salt and low-EDTA buffered solution. The preferred buffer is low TE (10 mM Tris, 0.1 mM EDTA, pH 8.0). The concentrations of Tris and EDTA must not exceed 10 mM and 0.1 mM, respectively. Avoid phosphate containing buffers.

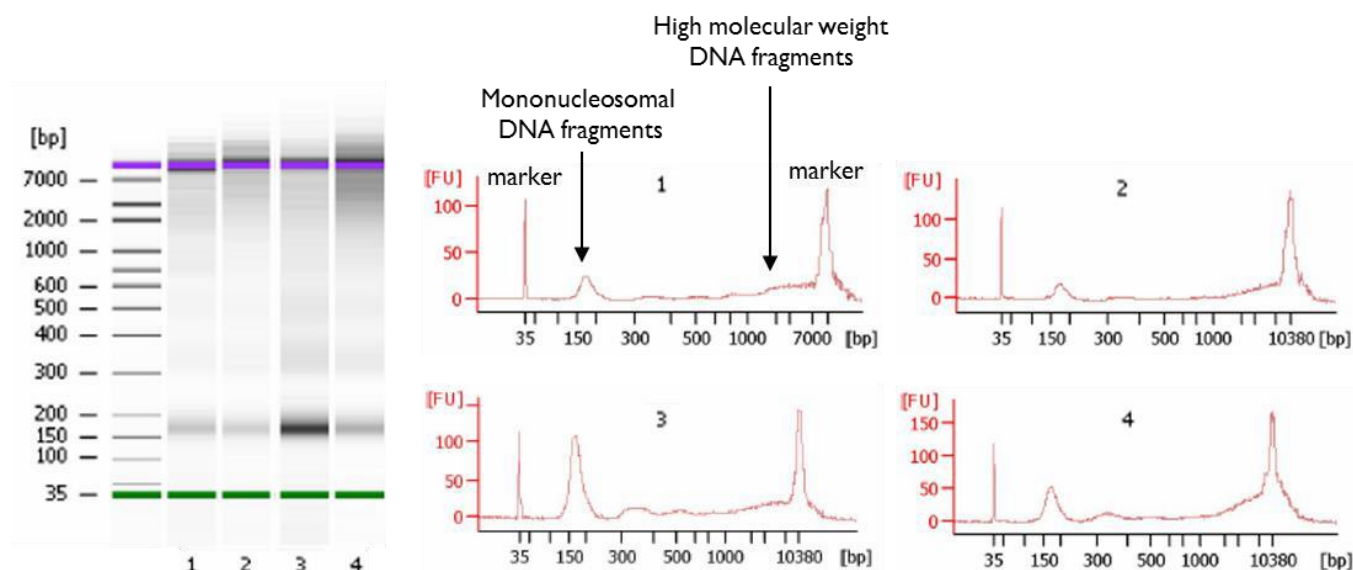
**6. Fragment Size**

The optimal DNA fragment size between 150 and 500 bp. The ThruPLEX DNA-Seq HV Kit is a ligation-based technology and adapters added during the process result in an approximately 140-bp increase in the size of each DNA template fragment. Library molecules with shorter inserts (200–300 bp) tend to cluster and amplify more efficiently on the Illumina flow cell. Depending on the application and requirements, the AMPure purification step following the final step (Library Amplification) can be replaced with a size-selection step to remove unwanted fragments.

**7. Using Cell-Free DNA from Plasma**

cfDNA isolated from plasma samples contains both high and low molecular weight DNA fragments (Figure 4). The composition and concentration of the isolated cfDNA differ from sample to sample and may vary depending on the isolation method used. The cfDNA species of most interest is the mononucleosomal DNA fragments of about 170 bp in length; the concentration can be determined using a fragment analyzer such as the Agilent Bioanalyzer.





**Figure 4. Fragment size distribution of cfDNA isolated from plasma.** cfDNA from four different human plasma samples was isolated and analyzed using the Agilent Bioanalyzer. Distinct features of the isolated cfDNA is a peak (or band) centered around 170 bp and higher molecular weight fragments.

## 8. Positive and Negative Controls

Include appropriate positive and negative controls in the experimental design to help verify that reactions proceed as expected. If the experimental samples contain any carryover contaminant(s) in the buffer, the downstream reactions may be impacted, and inclusion of controls would help elucidate such problems. Always prepare fresh dilutions of reference DNA (Control Fragmented Human gDNA, included in the kit). Include a negative control (No Template Control, NTC) with low TE buffer (10 mM Tris, 0.1 mM EDTA, pH 8.0) or nuclease-free water. The positive control and experimental samples should perform equivalently, while the NTC should not amplify.

## B. General Recommendations

### 1. Thermal Cycler Considerations

#### Thermal cycling and heated lid

Use a thermal cycler equipped with a heated lid that can handle 100- $\mu$ l reaction volumes. Set the temperature of the heated lid to 101°C–105°C to avoid sample evaporation during incubation and cycling.

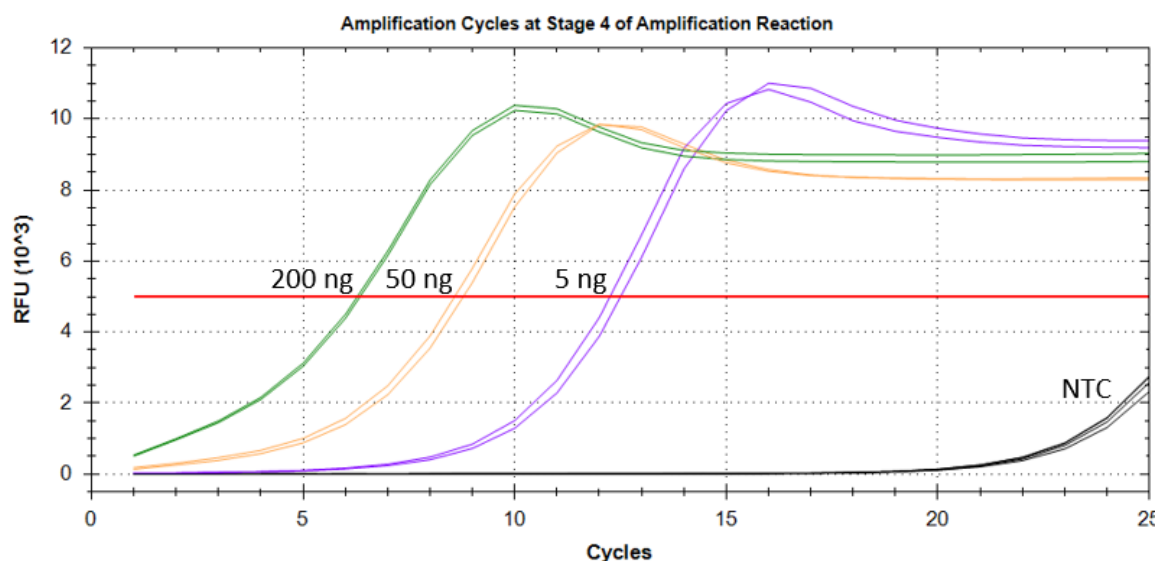
#### Thermal cycler ramp rates

We recommend a ramp rate of 3°C/s–5°C/s; higher ramp rates are not recommended and could impact the quality of the library.

#### Monitoring amplification during the Library Amplification Reaction

Amplification can be monitored using a real-time thermal cycler with the addition of fluorescent dyes (not provided with the kit, see Optional Materials in Section II.D) to the reaction (Figure 4). If a regular thermal cycler is used instead, there is no need to add the dyes; use an appropriate amount of nuclease-free water to prepare the Library Amplification Master Mix. In the absence of real-time monitoring, library amplification can be analyzed by gel or by analysis of an aliquot of the library using the Agilent Bioanalyzer (see Library Quantification, Section IV.B.3).

Depending on the real-time instrument used, select an appropriate calibration dye and mix with EvaGreen dye to prepare the dye mix (see Library Amplification Step, Section IV.A.3). For some real-time instruments, calibration dye may not be needed; please refer to the real-time thermal cycler instrument's manual.



**Figure 5. Example of real-time analysis of library amplification using ThruPLEX DNA-Seq HV.** A typical real-time amplification analysis of libraries prepared with ThruPLEX DNA-Seq HV Kit using 200 ng, 50 ng, or 5 ng of the provided positive control DNA compared to a No Template Control (NTC). Results were obtained using a CFX96 Touch Real-Time PCR Detection System (Bio-Rad Laboratories) with EvaGreen as the dye for detection and fluorescein as the calibration dye. The red line marks the midpoint of the linear phase of the amplification reaction curves and is used to determine the optimal number of amplification cycles at Stage 4 of the Library Amplification Reaction (Section IV.A). It is recommended to stay within one cycle above or below the optimal number of cycles. For example, for a 5-ng input, the optimal number of amplification cycles is  $12 \pm 1$  cycles or 11 to 13 cycles. The Relative Fluorescence Unit (RFU) values on the y-axis may vary based on the instrument used.

## 2. Preparation of Master Mixes

A master mix with appropriate buffers and enzymes must be prepared fresh at each workflow step, based on the number of reactions to be performed. Prepare ~10% excess of each master mix to allow for pipetting losses.

1. Transfer enzymes onto ice just prior to use and centrifuge briefly to collect contents at the bottom of the tube prior to use.
2. Thaw the buffers on ice, vortex briefly, and centrifuge prior to use.
3. Keep all components and master mixes on ice.
4. Once the master mix is prepared, thoroughly mix the contents several times with a pipette, while avoiding the introduction of air bubbles. Briefly centrifuge prior to dispensing into the PCR plate or tube(s).

The Library Synthesis Master Mix and Library Amplification Master Mix can be prepared during the last 15 min of the previous step's cycling protocol and kept on ice until use.

### 3. Indexing Reagents

We support use of the ThruPLEX HV index kits for use with ThruPLEX DNA-seq HV. The indexing modules are available to be purchased as a separate part or bundled with the core enzymatic components in a kit. They are available in 24-reaction or 96-reaction kit sizes and are available as unique dual indexes.

**NOTE:** The ThruPLEX DNA-Seq HV kit is **NOT** compatible with other versions of Takara Bio indexing kits, including those sold with previous versions of ThruPLEX kits. Contact Technical Support with any questions on compatibility.

Indexing Reagents consist of amplification primers containing Illumina-compatible indexes. Index sequences can be downloaded as .xlsx files at the ThruPLEX DNA-seq HV Product Page, under the Resources tab. Before starting the ThruPLEX DNA-seq HV Library Preparation Protocol (Section IV.A), refer to Appendix A for information on index sequences, Index Plate handling instructions, and multiplexing and index-pooling guidelines.

#### ThruPLEX HV UDI 24- and 96-reaction kits

Indexing Reagents are pre-dispensed and sealed in a linear barcoded Index Plate. The Index Plate is sealed with foil that can be pierced with a multichannel pipet tip to collect the required amount of index to assemble the reactions. Each well of the Index Plate contains sufficient volume for a single use. No more than four freeze/thaw cycles are recommended for the Index Plate.

### 4. Using Illumina Experiment Manager

Make sure the latest version of the Illumina Experiment Manager (IEM) is installed (version 1.18.1 or later). Prior to starting the ThruPLEX DNA-Seq HV Library Preparation Protocol (Section IV.A), create a Sample Sheet in the IEM to select and validate appropriate indexes to use in your experiments. Refer to Appendix A for guidelines on using the IEM to validate your index combinations.

### 5. Target Enrichment

ThruPLEX DNA-Seq HV is compatible with major target enrichment products. ThruPLEX DNA-Seq HV target enrichment protocols can be accessed through the Learning Center at [takarabio.com](http://takarabio.com).

## C. Safety Guidelines

Follow standard laboratory safety procedures and wear a suitable lab coat, protective goggles, and disposable gloves to ensure personal safety as well as to limit potential cross contaminations during the sample preparation and subsequent amplification reactions. For more information please refer to the appropriate Material Safety Data Sheets (MSDS) available online at [takarabio.com](http://takarabio.com).

## IV. Protocols

### A. Protocol: ThruPLEX DNA-Seq HV Library Preparation

#### 1. Template Preparation

**NOTE:** Assemble all reactions in thin-wall 96-well PCR plates or PCR tube(s) compatible with the thermal cycler and/or real-time thermal cycler used.

- Prepare samples as described below:
  - Samples:** Dispense 30 µl of fragmented doubled-stranded DNA, or cfDNA into each PCR tube or well of a PCR plate.
  - Positive control reactions using reference DNA:** If necessary, assemble reactions using 30 µl of the included Control Fragmented Human gDNA at an input amount comparable to the samples.
  - Negative control reactions/no-template controls (NTCs):** If necessary, assemble NTCs with 30 µl of nuclease-free water or TE buffer (e.g., 10 mM Tris, 0.1 mM EDTA, pH 8.0).

**NOTE:** The maximum volume of DNA cannot exceed 30 µl.

- Prepare **Template Preparation D Master Mix** as described in the table below for the desired number of reactions. Mix thoroughly with a pipette. Keep on ice until used:

| Template Preparation D Master Mix |           |                 |
|-----------------------------------|-----------|-----------------|
| Reagent                           | Cap color | Volume/reaction |
| PBD1                              | Blue      | 5 µl            |
| PED1                              | Blue      | 1 µl            |

**NOTE:** Prepare 10% excess to allow for pipetting losses.

- Assemble the **Template Preparation Reactions Mixture** as shown in the table below. To each 30 µl sample from Step 1 above, add 6 µl of the **Template Preparation Master Mix**.

| Template Preparation Reaction Mixture |                 |
|---------------------------------------|-----------------|
| Component                             | Volume/reaction |
| Sample or Control                     | 30 µl           |
| PBD1                                  | 5 µl            |
| PED1                                  | 1 µl            |
| <b>Total volume</b>                   | <b>36 µl</b>    |

- Mix thoroughly at least 10 times with a pipette set to 25 µl. Avoid introduction of air bubbles.
- Seal the PCR plate using an appropriate sealing film or tightly cap the tube(s).
- Centrifuge briefly to ensure the entire volume of the reaction is collected at the bottom of each well.
- Place the plate or tube(s) in a thermal cycler with heated lid set to 101°C–105°C. Perform the **Template Preparation Reaction** using the conditions in the table below:

| Template Preparation Reaction |                |
|-------------------------------|----------------|
| Temperature                   | Time           |
| 22°C                          | 25 min         |
| 55°C                          | 20 min         |
| 4°C                           | Hold for ≤2 hr |

## ThruPLEX® DNA-Seq HV User Manual

- After the thermal cycler reaches 4°C, remove the plate or tube(s) and centrifuge briefly before placing on ice.
- Proceed to the **Library Synthesis Step**.

**NOTE:** Following the Template Preparation step, spin down reaction and continue to Library Synthesis in the same plate or tube(s).

### 2. Library Synthesis

- Prepare **Library Synthesis D Master Mix** as described in the table below for the desired number of reactions. Mix thoroughly with a pipette. Keep on ice until used:

| Library Synthesis D Master Mix |           |                 |
|--------------------------------|-----------|-----------------|
| Reagent                        | Cap color | Volume/reaction |
| SBD1                           | White     | 1 µl            |
| SED1                           | White     | 4 µl            |

**NOTE:** Prepare 10% excess to allow for pipetting losses.

- Remove the seal on the plate or open the tube(s).
- Assemble the **Library Synthesis Reaction Mixture** as shown in the table below. To each well or tube, add 5 µl of the **Library Synthesis D Master Mix**.

| Library Synthesis Reaction Mix        |                 |
|---------------------------------------|-----------------|
| Component                             | Volume/reaction |
| Template Preparation Reaction Product | 36 µl           |
| SBD1                                  | 1 µl            |
| SED1                                  | 4 µl            |
| <b>Total volume</b>                   | <b>41 µl</b>    |

- Mix thoroughly at least 10 times with a pipette set to 25 µl.
- Seal the PCR plate using an appropriate sealing film or tightly cap the tube(s).
- Centrifuge briefly to collect the contents at the bottom of each well or tube.
- Return the plate or tube(s) to the thermal cycler with heated lid set to 101°C–105°C. Perform the **Library Synthesis Reaction** using the conditions in the table below:

| Library Synthesis Reaction |                  |
|----------------------------|------------------|
| Temperature                | Time             |
| 30°C                       | 40 min           |
| 4°C                        | Hold for ≤30 min |

- After the thermal cycler reaches 4°C, remove the plate or tube(s) and centrifuge briefly before placing on ice.
- Proceed to the **Library Amplification Step**.

**NOTE:** Following the Library Synthesis step, spin down reaction and continue to Library Amplification Reaction in the same plate or tube(s) maintained at 4°C.

**3. Library Amplification**

Multiple stages occur during the Library Amplification Reaction (see table in Step 8 below). During Stage 1 and Stage 2, the stem-loop adapters are extended and cleaved. **Proper programming of the thermal cycler is critical for these first two steps to be completed correctly, with no denaturation step occurring until Stage 3.** Illumina-compatible indexes are incorporated into the template library in Stage 4 and the resulting template is amplified; the number of cycles required at this stage is dependent on the amount of input DNA used. In stage 5 a final extension of the libraries occurs. Samples are cooled to 4°C in Stage 6, after which they are pooled and purified or stored at –20°C.

**Selection of the optimal number of cycles for library amplification (Stage 5):**

The number of PCR cycles required at Stage 5 of the Library Amplification Reaction is dependent on the amount of input DNA and thermal cycler used. Use the table below as a guide for selecting the number of PCR cycles.

| Stage 4 Amplification Guide |  |
|-----------------------------|--|
| Input DNA                   | Number of cycles required to generate a 500–1,000 ng library |
| 200 ng                      | 5–6  |
| 100 ng                      | 6–7  |
| 50 ng                       | 7–8  |
| 5 ng                        | 11–12  |

- **Optimization experiment:** Performing an optimization experiment to identify the appropriate number of PCR cycles needed is recommended. Use the desired amount of input DNA and allow the library amplification reaction to reach plateau. Determine the optimal number of amplification cycles by constructing PCR curves and identifying the midpoint of the linear phase as illustrated in Figure 4. Use the optimal amplification cycle number in the actual experiment for sequencing.
- **Yield:** The amount of amplified library can range from 100 ng to 1 µg depending upon many variables including sample type, fragmentation size, and thermal cycler used. When starting with Covaris-fragmented reference DNA with an average size of 200 bp and following the recommended number of amplification cycles, the typical yields range from 500 ng to 1000 ng.

**NOTE:** Over amplification could result in higher rate of PCR duplicates in the library.

**Reagents needed:**

| Reagent                     | Cap color   |
|-----------------------------|-------------|
| ABD1                        | Amber tube  |
| AED1                        | Violet      |
| Nuclease-Free Water         | Clear       |
| Fluorescent Dyes (optional) |             |
| Indexing Reagents           | Index Plate |

**NOTE:** It is critical to handle the Index Plate following the provided instructions to avoid cross contamination of indexes. If the entire Index Plate will not be used, please refer to Appendix A for Index Plate handling instructions. No more than four freeze/thaw cycles are recommended for the Index Plate.

## ThruPLEX® DNA-Seq HV User Manual

1. Prepare the Indexing Reagents described below:
  - Remove the Indexing Reagents from freezer and thaw for ten min on the bench top.
  - Spin the Indexing Reagents in a table top centrifuge to collect contents at the bottom of the well.
2. Prepare **Library Amplification D Master Mix** as described in the table below for the desired number of reactions. Mix thoroughly with a pipette. Keep on ice until used:

| Library Amplification D Master Mix           |            |                 |
|--|------------|-----------------|
| Reagent                                      | Cap color  | Volume/reaction |
| ABD1   | Amber tube | 42 µl           |
| AED1   | Violet     | 2 µl            |
| Nuclease Free Water (Plus Fluorescent Dyes*) | Clear      | 5 µl            |

### NOTES:

- Confirm ABD1 is fully thawed and thoroughly homogenized by heating briefly at 25°C and vortexing vigorously for 30 seconds.
- Prepare 10% excess to allow for pipetting losses.
- **If monitoring in real-time:** Fluorescence dyes\* (for detection and optical calibration) are added when monitoring amplification in real time during cycling. Please refer to the real-time PCR instrument's user manual for calibration dye recommendations. The volume of detection and calibration dyes plus nuclease-free water should not exceed 5 µl.
- **If not monitoring in real-time:** If a regular thermal cycler is used, there is no need to add the dyes; use 5 µl of nuclease-free water per reaction in the Library Amplification Master Mix.

3. Remove the seal on the PCR plate or open the tube(s).
4. Add 49 µl of the **Library Amplification D Master Mix** to each well or tube.
5. Add 10 µl of the appropriate Indexing Reagent to each well or tube:

### Index Plate Precautions:

- Make sure the two corner notches of the Index Plate are on the left, and the barcode label on the long side of the Index Plate is facing you.
- Thoroughly wipe the Index Plate seal with 70% ethanol and allow it to dry to prevent cross-contamination.

| Library Synthesis Reaction Mix               |                 |
|--|-----------------|
| Component                                    | Volume/reaction |
| Library Synthesis Reaction Product           | 41 µl           |
| ABD1   | 42 µl           |
| AED1   | 2 µl            |
| Nuclease Free Water (Plus Fluorescent Dyes*) | 5 µl            |
| ThruPLEX HV UDI                              | 10 µl           |
| <b>Total volume</b>                          | <b>100 µl</b>   |

6. Mix thoroughly a minimum of 10 times with a pipette set to 70 µl. Avoid introducing excessive air bubbles.
7. Seal the PCR plate using an appropriate sealing film or tightly cap the tube(s) and centrifuge briefly to collect the contents at the bottom of each well or tube.

**NOTE:** Use optical sealing film or caps if a real-time thermal cycler is used.

## ThruPLEX® DNA-Seq HV User Manual

- Return the plate or tube(s) to the thermal cycler with heated lid set to 101°C–105°C. Perform the **Library Amplification Reaction** using the conditions in the table below.

**CAUTION:** Ensure that the thermal cycler does not have a denaturing step programmed until Stage 3.

| Library Amplification Reaction |       |               |                  |  |
|--------------------------------|-------|---------------|------------------|--|
|                                | Stage | Temperature   | Time             | # Cycles                               |
| Extension & Cleavage           | 1     | 72°C          | 3 min            | 1                                      |
| Denaturation                   | 2     | 85°C          | 2 min            | 1                                      |
| Library Amplification          | 3     | 98°C          | 2 min            | 1                                      |
| Library Amplification          | 4     | 98°C<br>*68°C | 20 sec<br>75 sec | 5–12 (see Stage 4 Amplification Guide) |
| Final extension                | 5     | 68°C          | 5 min            | 1                                      |
| Hold                           | 6     | 4°C           | Hold             | 1                                      |

\*Acquire fluorescence data at this step, if monitoring amplification in real-time.

| Stage 4 Amplification Guide |  |
|-----------------------------|--|
| Input DNA                   | Number of cycles required to generate 500–1,000 ng library |
| 200 ng                      | 5–6  |
| 100 ng                      | 6–7  |
| 50 ng                       | 7–8  |
| 5 ng                        | 11–12  |

- Remove the PCR plate or tube(s) from the thermal cycler and centrifuge briefly to collect the contents to the bottom of each well.

**NOTE:** At this stage, samples can be processed for Next Generation Sequencing (NGS) immediately or stored frozen at –20°C for later processing. For instructions and recommendations on library pooling, purification, quantification, and sequencing, please refer to **Section IV.B**.

## B. Library Processing for Illumina Next-Generation Sequencing

### 1. Overview

This section contains guidelines for processing ThruPLEX DNA-Seq HV libraries for Illumina NGS. In some cases, recommended protocols are listed (Library Purification by AMPure XP beads) while in others, general guidelines are given. For more information, contact technical support at [technical\\_support@takarabio.com](mailto:technical_support@takarabio.com).

Libraries prepared from each sample will contain the specific indexes selected at the time of the amplification. needed. Once purified, the library should be quantified accurately prior to NGS to ensure efficient clustering on the Illumina flow cell. Instructions and recommendations on library purification and quantification, and quality are described in the following sections.



### 2. Library Purification by AMPure XP beads

AMPure XP is the recommended method of library purification. Do not use QIAquick cleanup or other silica-based filters for purification as this will result in incomplete removal of primers.

The ratio of AMPure XP beads to library DNA will determine the size-selection characteristics of the library. The ratio is also application dependent. For most NGS-based applications, a 1:1 bead to sample ratio is recommended. For more information, please refer to the vendor's recommendations on AMPure XP protocols for DNA purification.

#### Library purification reagents (supplied by the user)

| Reagent                              |
|--------------------------------------|
| AMPure XP beads                      |
| Magnetic rack for 200 µL strip tubes |
| Freshly prepared 80% (v/v) ethanol   |
| TE buffer, pH 8.0                    |

#### AMPure XP Protocol

##### NOTES:

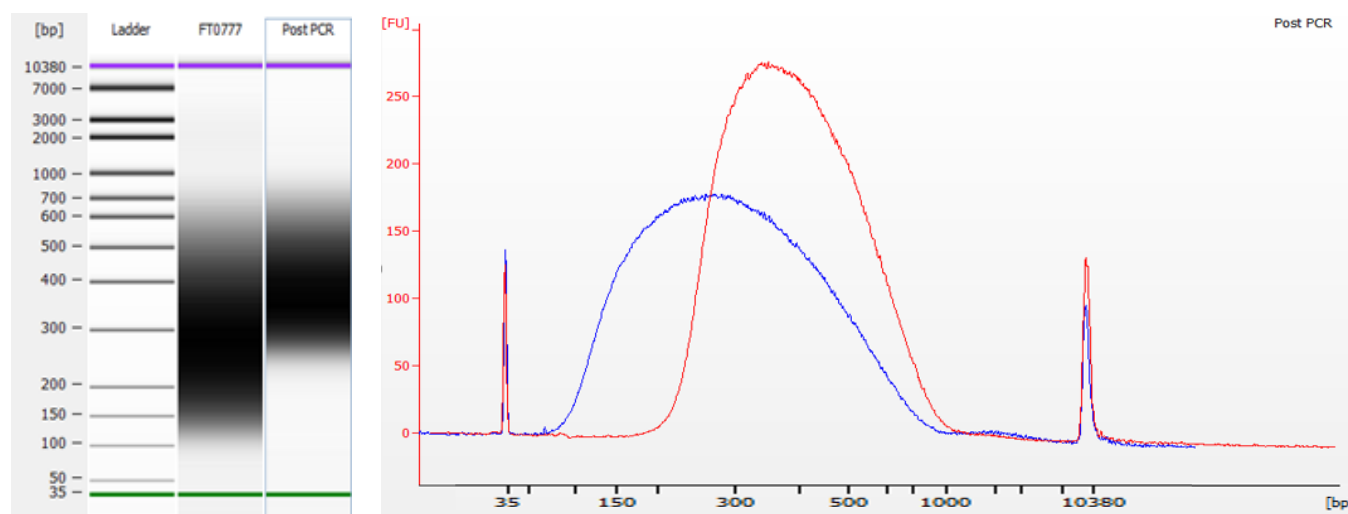
- It is important to bring all the samples and reagents to be used to room temperature.
  - Always use freshly prepared 80% (v/v) ethanol for Step 3 and Step 4 below.
  - Resuspend the AMPure XP reagent by gentle vortexing until no visible pellet is present at the bottom of the container.
1. In a 200 µl tube, mix 100 µl of AMPure XP reagent with 100-µl amplified library ensuring a 1:1 (v/v) ratio. Mix by pipette 10 times to achieve a homogeneous solution; incubate for 5 min at room temperature.
  2. Pulse-spin the sample(s) on a bench top centrifuge and place the tube in a magnetic stand. Wait for at least 2 min or until the beads are completely bound to the side of the tube(s) and the solution is clear.
  3. With the tube(s) in the magnetic stand and without disturbing the pellet, use a pipette to aspirate off and discard the supernatant. Add 200 µl of 80% (v/v) ethanol to the pellet and let stand for 30 seconds.
  4. With the tube(s) in the magnetic stand and without disturbing the pellet, use a pipette to aspirate off and discard the supernatant.
  5. Repeat Steps 3 and 4 for a total of two ethanol washes.
  6. Allow beads to air dry for no more than ~5 min—do not allow them to crack.
  7. Elute the DNA by re-suspending the beads with 50 µl of 1 x TE buffer, pH 8.0. Pulse-spin the sample(s) using a low speed, bench top centrifuge and place it into a magnetic stand and let the beads bind to the side of the tube(s) completely (for 2 min) until the solution is clear.
  8. While keeping the sample(s) in the magnetic stand, without disturbing the pellet, transfer the supernatant with a pipette into a new tube. If not used immediately, the purified library can be stored at -20°C.

## 3. Library Quantification and Quality Assessment

There are several approaches available for library quantification including real-time PCR with a library quantification kit for Illumina NGS libraries, such as Takara Library Quantification Kit (Takara Bio, Cat. No. 638324), fluorescence detection, using fluorescence detection-based methods, Qubit Fluorometer (Thermo Fisher Scientific), or Quant-iT PicoGreen dsDNA Assay Kit (Life Technologies), or using a fragment analyzer, such as the Agilent Bioanalyzer.

A fragment analyzer, such as the Agilent Bioanalyzer, can also be used to assess the quality of the libraries. We recommend diluting an aliquot of each library in TE buffer to ~5 ng/μL. Load a 1-μl aliquot of this diluted sample onto a Bioanalyzer high sensitivity DNA chip (Agilent Technologies, Cat. No. 5067-4626). Libraries prepared using the ThruPLEX DNA-Seq HV kit result in a size distribution of library fragments that is dependent on the input DNA fragment size (Figure 6).

**NOTE:** The adapters added during the ThruPLEX HV DNA-Seq library preparation process result in an approximately 140-base pair increase in the size of each library.



**Figure 6. Bioanalyzer analysis of libraries prepared using ThruPLEX DNA-Seq HV.** Libraries were prepared from 5-ng of the provided positive control DNA (Blue Trace) using the ThruPLEX DNA-Seq HV Kit. Post library amplification, libraries were purified following the AMPure XP protocol (IV.B.5). An aliquot of purified library was diluted to 5 ng/μl in TE buffer, and 1 μl of this diluted sample (Red Trace) was loaded onto a Bioanalyzer High Sensitivity DNA chip (Agilent Technologies).

## Appendix A. Indexing Reagents

### A. Overview

ThruPLEX DNA-Seq HV Kits are paired with ThruPLEX HV UDI kits containing unique dual-indexed PCR primers for amplification of indexed Illumina-compatible NGS libraries. These kits contain indexed PCR primers offering up to 96 unique dual indexes for multiplexing samples. The indexed PCR primers are supplied pre-dispensed in 96 well plates and are available in two formats: a set of 96 unique dual indexes (Cat. No. R400738) and a set of 24 unique dual indexes (Cat. No. R400739) that represents a subset of Cat. No. R400738. Each well of the dual index plate is for single use. All indexes have been functionally validated to work with Illumina sequencing systems using two- or four-channel chemistry for base calling. They have not been validated with systems using one-channel chemistry. Table 3 below summarizes the characteristics of the included Indexing Reagents, which consist of amplification primers containing Illumina-compatible indexes. Indexing Reagents should be stored at –20°C and should not be subjected to more than four freeze/thaw cycles.

**NOTE:** Indexing Reagents provided with ThruPLEX DNA-Seq HV kit cannot be substituted with indexing reagents from any other source.

### B. Components

Store all components at -20°C.

| Product Name          | Cat. No. | Concentration | Volume/tube |
|-----------------------|----------|---------------|-------------|
| ThruPLEX HV UDI 1-24* | R400739  | 12.5 µM       | 12 µl       |
| ThruPLEX HV UDI Set A | R400738  | 12.5 µM       | 12 µl       |

\*The indexes in the ThruPLEX HV UDI 1-24 kit are a subset of the ThruPLEX HV UDI Set A kit.

### C. ThruPLEX HV Unique Dual Index Sequences

The ThruPLEX HV unique dual indexes are 8-nt long and employ the “IDT for Illumina TruSeq® UD” i5 and i7 dual index sequences. An .xlsx file containing a full list of these indexes can be downloaded from our website.

Table V. ThruPLEX HV UDI - Set A plate layout.

|          | 1    | 2    | 3    | 4    | 5    | 6    | 7    | 8    | 9    | 10   | 11   | 12   |
|----------|------|------|------|------|------|------|------|------|------|------|------|------|
| <b>A</b> | U001 | U009 | U017 | U025 | U033 | U041 | U049 | U057 | U065 | U073 | U081 | U089 |
| <b>B</b> | U002 | U010 | U018 | U026 | U034 | U042 | U050 | U058 | U066 | U074 | U082 | U090 |
| <b>C</b> | U003 | U011 | U019 | U027 | U035 | U043 | U051 | U059 | U067 | U075 | U083 | U091 |
| <b>D</b> | U004 | U012 | U020 | U028 | U036 | U044 | U052 | U060 | U068 | U076 | U084 | U092 |
| <b>E</b> | U005 | U013 | U021 | U029 | U037 | U045 | U053 | U061 | U069 | U077 | U085 | U093 |
| <b>F</b> | U006 | U014 | U022 | U030 | U038 | U046 | U054 | U062 | U070 | U078 | U086 | U094 |
| <b>G</b> | U007 | U015 | U023 | U031 | U039 | U047 | U055 | U063 | U071 | U079 | U087 | U095 |
| <b>H</b> | U008 | U016 | U024 | U032 | U040 | U048 | U056 | U064 | U072 | U080 | U088 | U096 |

## ThruPLEX® DNA-Seq HV User Manual

Table VI. ThruPLEX HV UDI sequences.

| Index | i7 bases for sample sheet | i5 bases for sample sheet (MiSeq®, NovaSeq™, HiSeq® 2000/2500) | i5 bases for sample sheet (MiniSeq™, NextSeq®, HiSeq 3000/4000) |
|-------|---------------------------|--|---|
| U001  | CCGCGGTT                  | AGCGCTAG   | CTAGCGCT  |
| U002  | TTATAACC                  | GATATCGA   | TCGATATC  |
| U003  | GGACTTGG                  | CGCAGACG   | CGTCTGCG  |
| U004  | AAGTCCAA                  | TATGAGTA   | TACTCATA  |
| U005  | ATCCACTG                  | AGGTGCGT   | ACGCACCT  |
| U006  | GCTTGTC A                 | GAACATAC   | GTATGTTC  |
| U007  | CAAGCTAG                  | ACATAGCG   | CGCTATGT  |
| U008  | TGGATCGA                  | GTGCGATA   | TATCGCAC  |
| U009  | AGTTCAGG                  | CCAACAGA   | TCTGTTGG  |
| U010  | GACCTGAA                  | TTGGTGAG   | CTCACCAA  |
| U011  | TCTCTACT                  | CGCGGTTC   | GAACCGCG  |
| U012  | CTCTCGTC                  | TATAACCT   | AGGTTATA  |
| U013  | CCAAGTCT                  | AAGGATGA   | TCATCCTT  |
| U014  | TTGGACTC                  | GGAAGCAG   | CTGCTTCC  |
| U015  | GGCTTAAG                  | TCGTGACC   | GGTCACGA  |
| U016  | AATCCGGA                  | CTACAGTT   | AACTGTAG  |
| U017  | TAATACAG                  | ATATTACAC  | GTGAATAT  |
| U018  | CGGCGTGA                  | GCGCCTGT   | ACAGGCGC  |
| U019  | ATGTAAGT                  | ACTCTATG   | CATAGAGT  |
| U020  | GCACGGAC                  | GTCTCGCA   | TGCGAGAC  |
| U021  | GGTACCTT                  | AAGACGTC   | GACGTCTT  |
| U022  | AACGTTCC                  | GGAGTACT   | AGTACTCC  |
| U023  | GCAGAATT                  | ACCGGCCA   | TGGCCGGT  |
| U024  | ATGAGGCC                  | GTTAATTG   | CAATTAAC  |
| U025  | ACTAAGAT                  | AACCGCGG   | CCGCGGTT  |
| U026  | GTCGGAGC                  | GGTTATAA   | TTATAACC  |
| U027  | CTTGGTAT                  | CCAAGTCC   | GGACTTGG  |
| U028  | TCCAACGC                  | TTGGACTT   | AAGTCCAA  |
| U029  | CCGTGAAG                  | CAGTGGAT   | ATCCACTG  |
| U030  | TTACAGGA                  | TGACAAGC   | GCTTGTC A   |
| U031  | GGCATTCT                  | CTAGCTTG   | CAAGCTAG  |
| U032  | AATGCCTC                  | TCGATCCA   | TGGATCGA  |
| U033  | TACCGAGG                  | CCTGAACT   | AGTTCAGG  |
| U034  | CGTTAGAA                  | TTCAGGTC   | GACCTGAA  |
| U035  | AGCCTCAT                  | AGTAGAGA   | TCTCTACT  |
| U036  | GATTCTGC                  | GACGAGAG   | CTCTCGTC  |
| U037  | TCGTAGTG                  | AGACTTGG   | CCAAGTCT  |
| U038  | CTACGACA                  | GAGTCCAA   | TTGGACTC  |
| U039  | TAAGTGGT                  | CTTAAGCC   | GGCTTAAG  |
| U040  | CGGACAAC                  | TCCGGATT   | AATCCGGA  |
| U041  | ATATGGAT                  | CTGTATTA   | TAATACAG  |

## ThruPLEX® DNA-Seq HV User Manual

| Index | i7 bases for sample sheet | i5 bases for sample sheet (MiSeq, NovaSeq, HiSeq 2000/2500) | i5 bases for sample sheet (MiniSeq, NextSeq, HiSeq 3000/4000) |
|-------|---------------------------|---|---|
| U042  | GCGCAAGC                  | TCACGCCG  | CGGCGTGA  |
| U043  | AAGATACT                  | ACTTACAT  | ATGTAAGT  |
| U044  | GGAGCGTC                  | GTCCGTGC  | GCACGGAC  |
| U045  | ATGGCATG                  | AAGGTACC  | GGTACCTT  |
| U046  | GCAATGCA                  | GGAACGTT  | AACGTTCC  |
| U047  | GTTCCAAT                  | AATTCTGC  | GCAGAATT  |
| U048  | ACCTTGGC                  | GGCCTCAT  | ATGAGGCC  |
| U049  | ATATCTCG                  | ATCTTAGT  | ACTAAGAT  |
| U050  | GCGCTCTA                  | GCTCCGAC  | GTCGGAGC  |
| U051  | AACAGGTT                  | ATACCAAG  | CTTGGTAT  |
| U052  | GGTGAACC                  | GCGTTGGA  | TCCAACGC  |
| U053  | CAACAATG                  | CTTCACGG  | CCGTGAAG  |
| U054  | TGGTGGCA                  | TCCTGTAA  | TTACAGGA  |
| U055  | AGGCAGAG                  | AGAATGCC  | GGCATTCT  |
| U056  | GAATGAGA                  | GAGGCATT  | AATGCCTC  |
| U057  | TGCGGCGT                  | CCTCGGTA  | TACCGAGG  |
| U058  | CATAATAC                  | TTCTAACG  | CGTTAGAA  |
| U059  | GATCTATC                  | ATGAGGCT  | AGCCTCAT  |
| U060  | AGCTCGCT                  | GCAGAATC  | GATTCTGC  |
| U061  | CGGAACTG                  | CACTACGA  | TCGTAGTG  |
| U062  | TAAGGTCA                  | TGTGCTAG  | CTACGACA  |
| U063  | TTGCCTAG                  | ACCACTTA  | TAAGTGGT  |
| U064  | CCATTCGA                  | GTTGTCCG  | CGGACAAC  |
| U065  | ACACTAAG                  | ATCCATAT  | ATATGGAT  |
| U066  | GTGTCGGA                  | GCTTGCGC  | GCGCAAGC  |
| U067  | TTCCTGTT                  | AGTATCTT  | AAGATACT  |
| U068  | CCTTCACC                  | GACGCTCC  | GGAGCGTC  |
| U069  | GCCACAGG                  | CATGCCAT  | ATGGCATG  |
| U070  | ATTGTGAA                  | TGCATTGC  | GCAATGCA  |
| U071  | ACTCGTGT                  | ATTGGAAC  | GTTCCAAT  |
| U072  | GTCTACAC                  | GCCAAGGT  | ACCTTGGC  |
| U073  | CAATTAAC                  | CGAGATAT  | ATATCTCG  |
| U074  | TGGCCGGT                  | TAGAGCGC  | GCGCTCTA  |
| U075  | AGTACTCC                  | AACCTGTT  | AACAGGTT  |
| U076  | GACGTCTT                  | GGTTCACC  | GGTGAACC  |
| U077  | TGCGAGAC                  | CATTGTTG  | CAACAATG  |
| U078  | CATAGAGT                  | TGCCACCA  | TGGTGGCA  |
| U079  | ACAGGCGC                  | CTCTGCCT  | AGGCAGAG  |
| U080  | GTGAATAT                  | TCTCATTC  | GAATGAGA  |
| U081  | AACTGTAG                  | ACGCCGCA  | TGCGGCGT  |
| U082  | GGTCACGA                  | GTATTATG  | CATAATAC  |
| U083  | CTGCTTCC                  | GATAGATC  | GATCTATC  |

## ThruPLEX® DNA-Seq HV User Manual

| Index | i7 bases for sample sheet | i5 bases for sample sheet (MiSeq, NovaSeq, HiSeq 2000/2500) | i5 bases for sample sheet (MiniSeq, NextSeq, HiSeq 3000/4000) |
|-------|---------------------------|---|---|
| U084  | TCATCCTT                  | AGCGAGCT  | AGCTCGCT  |
| U085  | AGGTTATA                  | CAGTTCCG  | CGGAACTG  |
| U086  | GAACCGCG                  | TGACCTTA  | TAAGGTCA  |
| U087  | CTCACCAA                  | CTAGGCAA  | TTGCCTAG  |
| U088  | TCTGTTGG                  | TCGAATGG  | CCATTCTGA   |
| U089  | TATCGCAC                  | CTTAGTGT  | ACACTAAG  |
| U090  | CGCTATGT                  | TCCGACAC  | GTGTCCGA  |
| U091  | GTATGTTC                  | AACAGGAA  | TTCTGTGT  |
| U092  | ACGCACCT                  | GGTGAAGG  | CCTTCACC  |
| U093  | TACTCATA                  | CCTGTGGC  | GCCACAGG  |
| U094  | CGTCTGCG                  | TTCACAAT  | ATTGTGAA  |
| U095  | TCGATATC                  | ACACGAGT  | ACTCGTGT  |
| U096  | CTAGCGCT                  | GTGTAGAC  | GTCTACAC  |

## Appendix B. Troubleshooting Guide

Table VII. Troubleshooting Guide for the ThruPLEX Tag-Seq HV Kits.

| Problem   | Potential Cause   | Suggested Solutions   |
|---|---|---|
| Sample amplification curve looks like No Template Control (NTC) amplification curve or does not produce amplified product             | No input DNA added  | Quantitate input before using the kit   |
|   | Incorrect library template used (e.g., RNA, ssDNA)  | Adhere to DNA Sample Requirements (Section III.A)   |
| NTC amplification curve appears early or produces a yield similar to sample reaction products   | NTC contaminated with DNA   | Use a fresh control sample and check all reagents; replace kit if necessary.  |
|   |   | Clean area thoroughly and use PCR-dedicated plastics and pipettes.  |
| After purification of the amplified library, Bioanalyzer traces shows multiple peaks besides the markers                              | Input sample contains unevenly fragmented DNA of various sizes (e.g., plasma DNA)                               | If possible, quantify and check input DNA prior to using the kit.<br>Sequencing is still recommended.   |
| After purification of the amplified library, Bioanalyzer traces show broad peak(s) extending from 1,000 bp to greater than 10,000 bp. | Library was overamplified or the Bioanalyzer chip was overloaded. (This is common with high-sensitivity chips.) | Perform fewer PCR cycles during the Library Amplification Reaction.<br>For high-sensitivity chips, load ~1–5 ng/μl. Repeat the Bioanalyzer run. |

| Contact Us   |  |
|--|--|
| Customer Service/Ordering  | Technical Support  |
| tel: 800.662.2566 (toll-free)  | tel: 800.662.2566 (toll-free)  |
| fax: 800.424.1350 (toll-free)  | fax: 800.424.1350 (toll-free)  |
| web: <a href="http://takarabio.com">takarabio.com</a>                      | web: <a href="http://takarabio.com">takarabio.com</a>  |
| e-mail: <a href="mailto:ordersUS@takarabio.com">ordersUS@takarabio.com</a> | e-mail: <a href="mailto:technical_support@takarabio.com">technical_support@takarabio.com</a> |

## Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Takara Bio USA, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at [takarabio.com](http://takarabio.com). It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

**©2020 Takara Bio Inc. All Rights Reserved.**

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at [takarabio.com](http://takarabio.com).

This document has been reviewed and approved by the Quality Department.