

NEBNext[®] Immune Sequencing Kit (Human)

NEB #E6320S/L

24/96 reactions

Version 2.0_12/21

Table of Contents

Overview	2
Applications.....	5
Protocol.....	6
Kit Components.....	15
Checklist.....	17
Revision History	20

The Kit Includes

NEB #E6320S contains E2628S (Pouch with BCR and TCR primers), #E6325-1 (Box 1 of 2), and #E6325-2 (Box 2 of 2) and is sufficient for preparation of up to 24 reactions. NEB #E6320L contains E2628L (Pouch with BCR and TCR primers), #E6325-3 (Box 1 of 2), and #E6325-4 (Box 2 of 2) and is sufficient for preparation of up to 96 reactions. All reagents should be stored at -20°C , 4°C or room temperature as indicated below. Colored bullets represent the color of the cap of the tube containing the reagent.

NEB #E2628S/L, NEBNext Immune Sequencing Primers (Human)

Pouch: Store at -20°C .

- (blue) NEBNext IS BCR Primers (Human)
- (blue) NEBNext IS TCR Primers (Human)

NEB #E6325-1, NEBNext Immune Sequencing Kit RT and PCR Reagents (NEB #E6320S, 24 reactions)

Box 1 of 2: Store at -20°C .

- (lilac) NEBNext IS RT Buffer (4X)
- (lilac) NEBNext IS RT Primer
- (lilac) NEBNext IS TS Oligos
- (lilac) NEBNext First Strand Synthesis Enzyme Mix
- (lilac) dNTP Solution Mix
- (lilac) NEBNext Cell Lysis Buffer
- (blue) Q5[®] Reaction Buffer (5X)
- (blue) Q5 Hot Start High-Fidelity DNA Polymerase
- (blue) NEBNext i701-i706 Primers
- (blue) NEBNext i501-i504 Primers
- (blue) NEBNext IS PCR2 Universal Primer
- (white) Nuclease-free Water
- (white) TE (0.1X)

NEB #E6325-3, NEBNext Immune Sequencing Kit RT and PCR Reagents (NEB #E6320L, 96 reactions)

Box 1 of 2: Store at –20°C.

- (lilac) NEBNext IS RT Buffer (4X)
- (lilac) NEBNext IS RT Primer
- (lilac) NEBNext IS TS Oligos
- (lilac) NEBNext First Strand Synthesis Enzyme Mix
- (lilac) dNTP Solution Mix
- (lilac) NEBNext Cell Lysis Buffer
- (blue) Q5 Reaction Buffer (5X)
- (blue) Q5 Hot Start High-Fidelity DNA Polymerase
- (blue) NEBNext i701-i712 Primers
- (blue) NEBNext i501-i508 Primers
- (blue) NEBNext IS PCR2 Universal Primer
- (white) Nuclease-free Water
- (white) TE (0.1X)

NEB #E6325-2/#E6325-4, NEBNext Immune Sequencing Kit cDNA Purification Reagents

Box 2 of 2: Store at 4°C. Do not freeze.

- (white) Streptavidin Magnetic Beads
- (white) NEBNext IS Bind and Wash Buffer (2X)
- (white) 0.1% Tween
- (white) NEBNext Sample Purification Beads*

*This component should be stored at room temperature.

Required Materials Not Included

- SYBR® Green (Life Technologies #S7563)
- 80% Ethanol (freshly prepared)
- DNA LoBind® Tubes (Eppendorf #022431021)
- Magnetic rack/stand (NEB #S1515, Alpaqua®, cat. #A001322 or equivalent)
- Thermal cycler
- qPCR machine and associated 96 well plates
- Bioanalyzer®/TapeStation® and associated consumables
- MiSeq® and v3 600 cycle sequencing reagent
- Nuclease-free Water
- Tube rotating mixer
- Magnet stand or equivalent (for washing beads only) (NEB #S1506)
- DNase RNase-free PCR strip tubes (USA Scientific 1402-1708)

Overview

Immune repertoire sequencing is a powerful tool that can be used to analyze the immune response to diseases and pathogens from current and past exposures. Areas of particular interest include characterization of autoimmune diseases, oncology, discovery of neutralizing antibodies against infectious disease, or use as a diagnostic tool for residual disease detection.

In order to create the incredible genetic diversity required for antigen recognition by B and T Cells, lymphocytes construct unique antigen receptors via a process known as V(D)J recombination. In this process, each cell selects a single V, D, and J gene segment through genetic recombination, introducing additional non-germline-encoded nucleotides at the junctions. This process generates immune receptor diversity, the majority of which is encoded in the heavy chain complementarity determining region 3 (CDR3) (Figure 1).

Recent improvements in the read length and throughput of next-generation sequencing (NGS) platforms have resulted in a rise in the popularity of immune repertoire sequencing. The complex structure and sequence diversity of antibody-encoding genes have provided significant challenges to the development of a simple and reliable method to sequence immune receptor sequences.

The NEBNext Immune Sequencing Library Preparation Kit (Figure 2) provides all the components required for enrichment and sequencing of the B Cell Receptor (BCR) and T Cell Receptor (TCR) RNA transcripts present in a sample. NEBNext Immune Sequencing Kit has been developed and optimized to provide accurate sequencing of full-length immune transcript repertoires of B-cells and T-cells. This allows for exhaustive somatic mutation profiling across complete V, D and J segments, full isotype information analysis (IgM, IgD, IgG, IgA and IgE), and also allows for synthesis and expression of complete antibody chains for downstream immunological assays. In addition, this approach allows characterization of TRA and TRB chains. The method uses unique molecular identifier scheme specifically designed to barcode each mRNA molecule, allowing PCR copies derived from an individual mRNA to be collapsed into a single consensus sequence. This improves sequence accuracy by resolving PCR and sequencing errors. This also eliminates PCR bias, thus allowing for quantitative digital molecule counting.

Defining features of the Immune Sequencing method include:

- a) Full-length variable sequences are generated (including isotype information), allowing downstream antibody synthesis and functional characterization not possible with approaches sequencing only the CDR3 region.
- b) Variable region primers are not required, which reduces primer pool complexity and allows for the unbiased and simultaneous recovery of B Cell and T Cell receptor transcripts.
- c) Unique molecular barcoding approach minimizes PCR biases and improves sequencing accuracy by allowing a consensus to be generated from duplicate sequencing reads originating from the same transcript. Moreover, unique barcoding enables accurate quantitation of each clone present in the sample.
- d) High target capture efficiency allows for immune repertoire sequencing and analysis from sub-microgram quantities of total RNA.

Figure 1. Simplified representation of the structure of BCR or TCR showing the outcome of V(D)J recombination in mature lymphocytes.

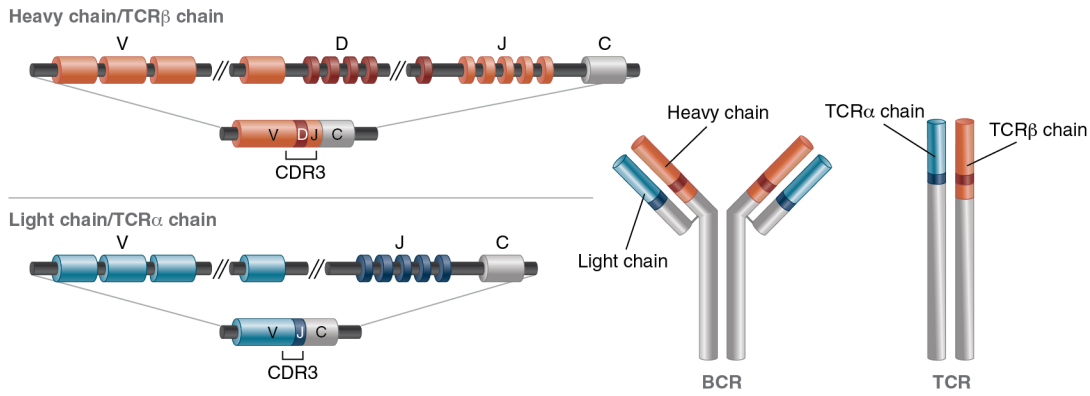
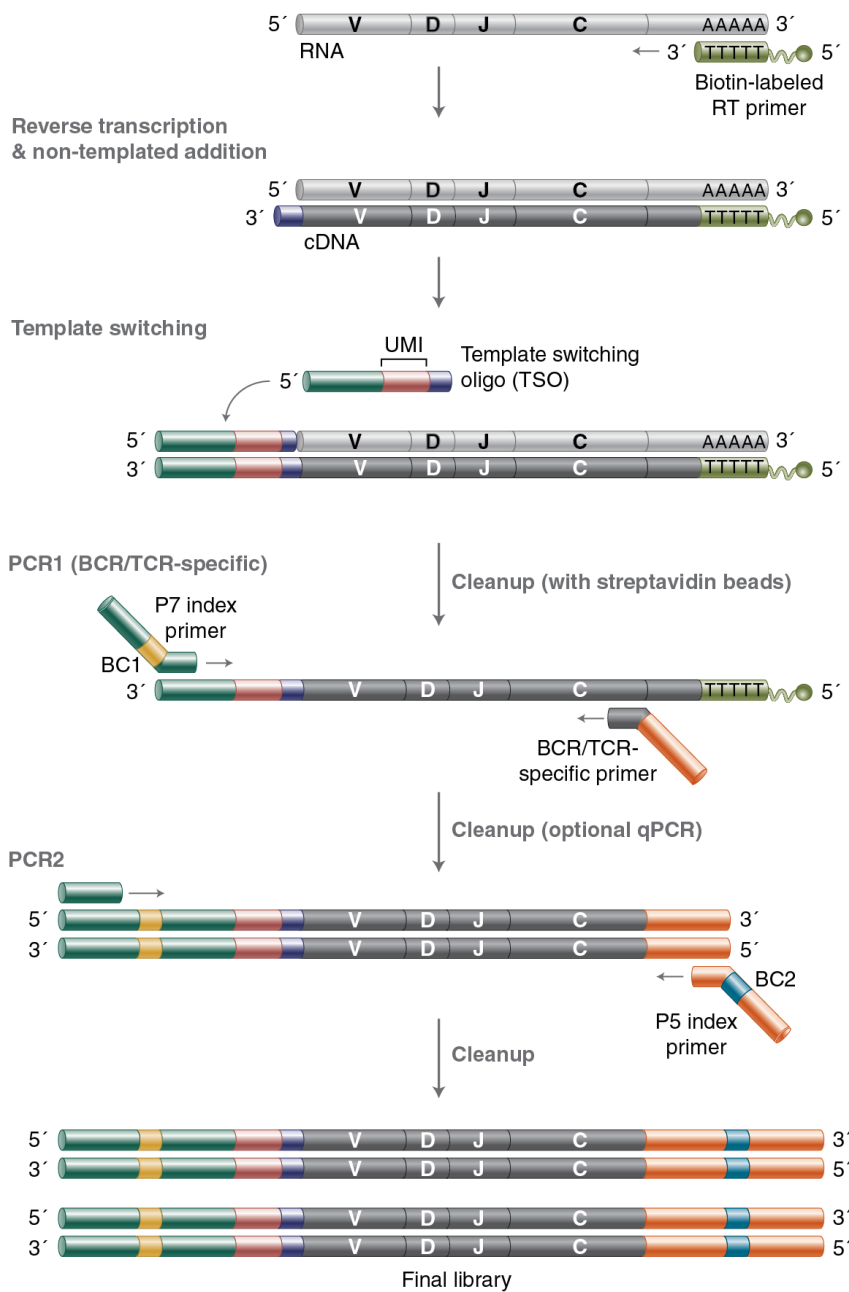


Figure 2. NEBNext Immune Sequencing Library Preparation Kit Workflow



Step I—Prepare target RNA

Immune Sequencing does not require mRNA isolation prior to library construction. Total RNA can be prepared from any sample or tissue known to contain B Cells or T Cells, such as peripheral blood mononuclear cells, bone marrow, lymph nodes, or any other tissue in which infiltrating lymphocytes are present. Cells or tissues can be frozen, but intact, high quality RNA is required for optimal performance.

Step II—Perform molecular-tagging reverse transcription

During this 50-minute reaction full-length mRNA is reverse transcribed and the 3' end of each cDNA is tagged with a DNA adaptor containing a unique molecular identifier (UMI) barcode.

Step III—Purify with Streptavidin Magnetic Beads

This 30-minute cleanup step removes all salts, reagents and unused cDNA adaptor oligonucleotides (TSO and RT Primers) allowing clean downstream amplification steps.

Step IV—Primary amplification of target fragments

This primary PCR step uses one primer which binds the universal cDNA adaptor and a primer mix targeting constant regions of the desired immune receptors. Products are then purified with Sample Purification Beads.

Step V—Secondary PCR cycle optimization with real-time PCR

This optional real-time PCR allows accurate assessment of target quantity leading to an optimized choice of PCR cycles for the final library PCR.

Step VI—Secondary PCR including Illumina® adaptor addition

This secondary PCR adds additional Illumina adaptor sequences to the target fragments (including Illumina multiplexing index sequences if desired). After Sample Purification Bead cleanup the libraries are ready for Bioanalyzer or TapeStation analysis, quantification and sequencing on the MiSeq platform.

Applications

The NEBNext Immune Sequencing Kit (Human) contains enzymes and buffers that are ideal to convert a small amount of human RNA input into indexed libraries for next-generation sequencing on the Illumina® MiSeq platform (Illumina, Inc). The workflow of NEBNext Immune Sequencing (Human) Kit is very user-friendly and fast with minimal hands-on time. Each of these components must pass rigorous quality control standards and are lot controlled, both individually and as a set of reagents.

For larger volume requirements, customized and bulk packaging is available by purchasing through the OEM/Bulks department at NEB. Please contact OEM@neb.com for further information.

Protocol for NEBNext Immune Sequencing Kit (Human)

Symbols



This is a point where you can safely stop the protocol and store the samples prior to proceeding to the next step in the protocol.



This caution sign signifies a step in the protocol that has two paths leading to the same end point.



Colored bullets indicate the cap color of the reagent to be added.

RNA Sample Recommendations: The RNA sample should be free of salts (e.g., Mg^{2+} , or guanidine salts) or organics (e.g., phenol and ethanol). High quality RNA with a RIN > 7 is recommended for optimal performance.

Starting Material: 10 ng–1 μ g enriched B cell or T Cell or B cell and T Cell total RNA, or 10 ng–1 μ g total RNA from peripheral blood mononuclear cells (PBMCs). Note: The maximum volume of RNA input is 9 μ l.

1. NEBNext Immune Sequencing Reverse Transcription and cDNA Synthesis

1.1. Mix the following components in a sterile nuclease-free tube:

COMPONENT	VOLUME
• (lilac) NEBNext IS RT Buffer (4X)	5 μ l
• (lilac) NEBNext IS RT Primer	1 μ l
• (lilac) NEBNext IS TS Oligos	1 μ l
• (lilac) dNTP Solution Mix	2 μ l
• (lilac) NEBNext First Strand Synthesis Enzyme Mix	2 μ l
• (lilac) NEBNext Cell Lysis Buffer	0.5 μ l
Total RNA	1-9 μ l
○ (white) Nuclease-free Water	Variable
Total Volume	20.5 μ l

1.2. Set a 100 μ l or 20 μ l pipette to 15 μ l and then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube. Note: It is important to mix well. The presence of a small amount of bubbles will not interfere with the performance.

1.3. Place in a thermal cycler, with the heated lid set to $\geq 80^{\circ}C$, and run the following program:

CYCLE STEP	TEMP	TIME
Incubation	42 $^{\circ}C$	40 minutes
Inactivation	70 $^{\circ}C$	10 minutes
Hold	4 $^{\circ}C$	∞

2. Purify the cDNA with Streptavidin Magnetic Beads

2.1. Prepare Bind and Wash Buffer (1X) to be used in Step 2.11.

- Dilute NEBNext IS Bind and Wash Buffer (2X) to 1X with an equal volume of nuclease-free water (not provided).
- For each sample, you will need 100 μ l Bind and Wash Buffer (1X).

2.2. Vortex Streptavidin Magnetic Beads to resuspend. Aliquot the total amount of Streptavidin Magnetic Beads needed (15 μ l of Streptavidin Magnetic Beads per sample) into a clean RNase-free 1.5 ml tube.

2.3. Place the tube on a magnetic rack at room temperature. Once the solution is clear (~2 minutes) carefully remove and discard the supernatant without disturbing the bead pellet.

2.4. Remove the tube from the magnet and wash the beads by adding 200 μ l of undiluted NEBNext IS Bind and Wash Buffer (2X). Vortex briefly to resuspend the beads and quickly spin the tube in a microcentrifuge to collect any sample on the sides of the tube.

2.5. Place the tube on a magnetic rack. Once the solution is clear, carefully remove and discard the supernatant without disturbing the bead pellet.

- 2.6. Repeat Steps 2.4–2.5 once for a total of two washes.
- 2.7. Remove the tube from the magnetic rack. Resuspend the beads in 20 µl of undiluted NEBNext IS Bind and Wash Buffer (2X) for each sample (e.g., for 4 samples, add 4 x 20 = 80 µl NEBNext IS Bind and Wash Buffer (2X). Vortex briefly to resuspend the beads.
- 2.8. Add 20 µl beads to the cDNA products (20.5 µl) from Step 1.3. Vortex briefly to mix.
- 2.9. Place the tube on a rotator for 15 minutes at room temperature.
- 2.10. Quickly spin down the tube to collect any sample on the sides of the tube and place the tube on a magnetic rack. Once the solution is clear, carefully remove and discard the supernatant without disturbing the bead pellet.
- 2.11. Remove the tube from the magnetic rack. Add 100 µl diluted Bind and Wash Buffer (1X) (from Step 2.1). Vortex briefly to resuspend the beads (Note: vortexing is better than flicking the tube, as it minimizes foaming).
- 2.12. Quickly spin down the tube to collect any sample on the sides and place the tube on a magnetic rack. Once the solution is clear, carefully remove and discard the supernatant without disturbing the bead pellet.
- 2.13. Remove the tube from the magnetic rack and add 100 µl of 0.1% Tween. Vortex briefly to resuspend the beads (Note: This step is critical to remove all the salt from the beads before elution).
- 2.14. Quickly spin down the tube to collect any sample on the sides of the tube and place the tubes on a magnetic rack. Once the solution is clear, carefully remove and discard the supernatant without disturbing the bead pellet. With a 20 µl pipette tip, carefully remove any residual liquid left at the bottom of the tube without disturbing the beads.
- 2.15. Remove the tube from the magnet and add 23 µl of 0.1% Tween. Vortex briefly to resuspend the beads and quickly spin in a microcentrifuge to collect any sample on the sides of the tube.
- 2.16. Place the tube in a thermal cycler, with the heated lid set at 105°C, and run the following program:

CYCLE STEP	TEMP	TIME
Incubation	95°C	3 minutes
Hold	25°C	∞

- 2.17. Remove the tube from the thermal cycler, vortex briefly, spin down the contents and place on the magnetic rack.
- 2.18. Once the solution is clear, transfer 21 µl of the supernatant into a new tube and discard the beads.



Safe Stopping Point: It is safe to store the cDNA at -20°C overnight.

3. PCR1 for VDJ Region Amplification and Index Barcode Incorporation



Use Option 3.1A for enriching B Cell Receptor (BCR) chains.

Use Option 3.1B for enriching T Cell Receptor (TCR) chains.

Use Option 3.1C for enriching both BCR chains and TCR chains.

- 3.1. Add the following components to a sterile strip tube:

3.1A. PCR1 for Enriching B Cell Receptor Chains

COMPONENT	VOLUME
Purified cDNA (Step 2.18)	21 µl
• (blue) Q5 Reaction Buffer (5X)	10 µl
• (lilac) dNTP Solution Mix	1 µl
• (blue) Q5 Hot Start High-Fidelity DNA Polymerase	1 µl
• (blue) NEBNext IS BCR Primers (Human)	2 µl
• (blue) NEBNext i7 Index Primer*	1 µl
○ (white) Nuclease-free Water	14 µl
Total Volume	50 µl

*Use only one i7 Primer per sample

3.1B. PCR1 for Enriching T Cell Receptor Chains

COMPONENT	VOLUME
Purified cDNA (Step 2.18)	21 µl
• (blue) Q5 Reaction Buffer (5X)	10 µl
• (lilac) dNTP Solution Mix	1 µl
• (blue) Q5 Hot Start High-Fidelity DNA Polymerase	1 µl
• (blue) NEBNext IS TCR Primers (Human)	2 µl
• (blue) NEBNext i7 Primer*	1 µl
○ (white) Nuclease-free Water	14 µl
Total Volume	50 µl

*Use only one i7 index primer per sample

3.1C. PCR1 for Enriching Both B Cell Receptor Chains and T Cell Receptor Chains

COMPONENT	VOLUME
Purified cDNA (Step 2.18)	21 µl
• (blue) Q5 Reaction Buffer (5X)	10 µl
• (lilac) dNTP Solution Mix	1 µl
• (blue) Q5 Hot Start High-Fidelity DNA Polymerase	1 µl
• (blue) NEBNext IS BCR Primers (Human)	2 µl
• (blue) NEBNext IS TCR Primers (Human)	2 µl
• (blue) NEBNext i7 Index Primer*	1 µl
○ (white) Nuclease-free Water	12 µl
Total Volume	50 µl

*Use only one i7 Primer per sample

- 3.2. Set a 100 µl or 200 µl pipette to 40 µl and then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube.
- 3.3. Place the tube on a thermal cycler and perform PCR amplification using the following PCR cycling program:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	60 seconds	1
Denaturation	98°C	10 seconds	12
Annealing	64°C	30 seconds	
Extension	72°C	30 seconds	
Final Extension	72°C	60 seconds	1
Hold	4°C	∞	

4. Cleanup of PCR1 Amplification

- 4.1. Vortex NEBNext Sample Purification Beads to resuspend. SPRIselect and AMPure® XP Beads can be used as well. If using AMPure XP Beads, please allow the beads to warm to room temperature for at least 30 minutes before use.
- 4.2. Add 50 µl of resuspended NEBNext Sample Purification Beads to each 50 µl PCR reactions. Mix well by pipetting or brief vortexing.
- 4.3. Incubate for 5 minutes at room temperature.
- 4.4. Quickly spin the tube in a microcentrifuge and place it on a magnetic rack. Once the solution is clear (~5 minutes) carefully remove and discard the supernatant without disturbing the bead pellet. (**Caution: do not discard beads**).

- 4.5. Wash beads by adding 200 µl of freshly prepared 80% ethanol to the tube while in the magnetic rack. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant.
- 4.6. Repeat Step 4.5 once for a total of two wash steps. After supernatant is discarded, use a 20 µl pipette tip to carefully remove any residual liquid left at the bottom of the tube without disturbing the beads.
- 4.7. Air dry the beads for up to 5 minutes while the tube is on the magnetic stand with the lid open.
- Caution: Do not over dry the beads. This may result in lower recovery of DNA target.**
- 4.8. Elute DNA target from beads by adding 25 µl TE (0.1X) to beads. Mix well on a vortex mixer or by pipetting up and down at least 10 times. Incubate for at least 2 minutes at room temperature. Quickly spin the tube and place it on a magnetic rack to separate beads from supernatant.
- 4.9. After the solution is clear (about 2 minutes), carefully transfer 23 µl supernatant to a new PCR tube.



Safe Stopping Point: It is safe to store the PCR product overnight at -20°C.

5. qPCR for PCR2 Cycle Optimization

The qPCR step is optional; it is only used to identify the optimal final PCR cycling conditions, as each sample can contain varying amounts of amplifiable molecules. One can achieve a similar result by manually taking an aliquot of the PCR reaction every 5 cycles and visualizing each aliquot on Bioanalyzer or TapeStation to identify the optimal cycling number that yields the highest product with the lowest background. The number of cycles should be high enough to provide sufficient library fragments for a successful sequencing run, but low enough to avoid PCR artifacts and over-cycling (high molecular weight fragments on Bioanalyzer). Typically, most samples will amplify best in the range of 6–18 cycles.

- 5.1. Add the following components to a sterile strip tube:

COMPONENT	VOLUME
Purified PCR1 DNA (Step 4.9)	5 µl
• (blue) Q5 Reaction Buffer (5X)	5 µl
• (lilac) dNTP Solution Mix	0.5 µl
• (blue) Q5 Hot Start High-Fidelity DNA Polymerase	0.5 µl
• (blue) NEBNext IS PCR2 Universal Primer	1 µl
• (blue) NEBNext i5 Primer*	1 µl
SYBR Green (20X)**	0.25 µl
○ (white) Nuclease-free Water	11.75 µl
Total Volume	25 µl

* Use one i5 Primer per PCR reaction

** 20X working stock of SYBR Green can be made by diluting 1 µl (10,000X) SYBR Green (Life Technologies #S7563) into 399 µl DMSO and stored at 4°C for up to a month. You only need 0.2 x SYBR final in the qPCR reaction.

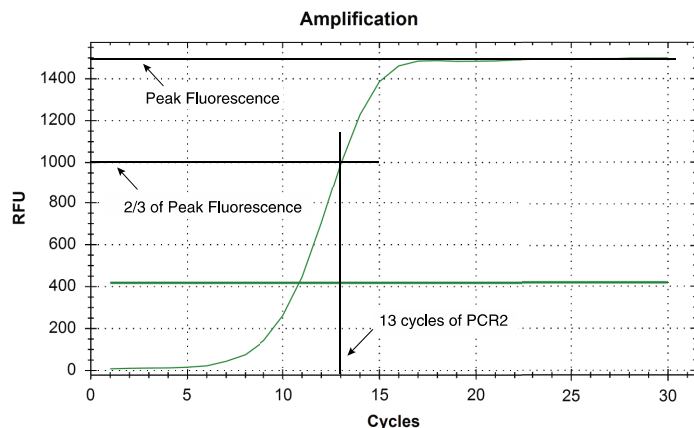
- 5.2. Set a 100 µl or 20 µl pipette to 20 µl and then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube.
- 5.3. Place the tube on a qPCR instrument and run the following PCR cycling program:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	60 seconds	1
Denaturation	98°C	10 seconds	30
Annealing and Extension	72°C	30 seconds (fluorescence on*)	
Hold	4°C	∞	

* Select the SYBR channel.

- 5.4. The amplification curve for each sample is used to determine the optimal number of PCR cycles to use during PCR2 (Section 6). The PCR cycle number is approximated by determining the cycle number (round down to the nearest whole number) corresponding to 2/3 of the peak fluorescence (see example in Figure 3; in this example 13 PCR cycles will be used in Section 6).

Figure 3. Determining cycle number for PCR2 using amplification curve of sample from a qPCR run.



6. PCR2 for Final Library

- 6.1. Add the following components to a sterile strip tube:

COMPONENT	VOLUME
Purified PCR1 DNA (Step 4.9)	10 μ l
• (blue) Q5 Reaction Buffer (5X)	10 μ l
• (lilac) dNTP Solution Mix	1 μ l
• (blue) Q5 Hot Start High-Fidelity DNA Polymerase	1 μ l
• (blue) NEBNext IS PCR2 Universal Primer	2 μ l
• (blue) NEBNext i5 Index Primer*	2 μ l
○ (white) Nuclease-free Water	24 μ l
Total Volume	50 μ l

* Use one i5 Primer per PCR reaction

- 6.2. Set a 100 μ l pipette to 40 μ l then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube.
- 6.3. Place the tube on a thermal cycler and run the following PCR cycling program:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	60 seconds	1
Denaturation	98°C	10 seconds	X*
Annealing and Extension	72°C	30 seconds	
Hold	4°C	∞	

* Use number of cycles determined from Step 5.4. If qPCR was not done, see Table 1 for recommended number of PCR cycles. In this case the number of PCR cycles should be chosen based on input amount and sample type. The number of cycles should be high enough to provide sufficient library fragments for a successful sequencing run, but low enough to avoid PCR artifacts and over-cycling (high molecular weight fragments on Bioanalyzer).

Table 1.

For enriched B Cell or T Cell RNA input, PCR2 cycle number is recommended as in the table:

TOTAL RNA INPUT	PCR1 ASSAY	PCR2 CYCLE NUMBER
B Cell RNA 1,000 ng	BCR	6-7
B Cell RNA 100 ng	BCR	9-10
B Cell RNA 10 ng	BCR	13-14
T Cell RNA 1,000 ng	TCR	10-11
T Cell RNA 100 ng	TCR	12-13
T Cell RNA 10 ng	TCR	17-18

Note: if B and T cells are combined, choose a PCR2 cycle number in between the recommendations for B cell and T cell RNA.

7. Cleanup of PCR2 Amplification to Obtain Final Libraries

- 7.1. Vortex NEBNext Sample Purification Beads to resuspend. SPRIselect and AMPure XP Beads can be used as well. If using AMPure XP Beads, please allow the beads to warm to room temperature for at least 30 minutes before use.
- 7.2. Add 50 μ l of TE (0.1X) to each 50 μ l PCR reaction to obtain 100 μ l mixture, and then add 50 μ l resuspended NEBNext Sample Purification Beads to the 100 μ l mixture. Mix well by pipetting or brief vortexing.
- 7.3. Incubate for 5 minutes at room temperature.
- 7.4. Quickly spin the tube in a microcentrifuge and place it on a magnetic rack. Once the solution is clear (~5 minutes) carefully transfer the 150 μ l supernatant to a new tube without disturbing the bead pellet. **Save the supernatant and discard beads.**
- 7.5. Add 15 μ l of resuspended NEBNext Sample Purification Beads to supernatant. Mix well by pipetting or brief vortexing.
- 7.6. Incubate for 5 minutes at room temperature.
- 7.7. Quickly spin the tube in a microcentrifuge and place it on a magnetic rack. Once the solution is clear (~5 minutes) carefully remove and discard the supernatant without disturbing the bead pellet. **Do not discard beads.**
- 7.8. Wash beads by adding 200 μ l of freshly prepared 80% ethanol to the tube while in the magnetic rack. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant.
- 7.9. Repeat Step 7.8 once for a total of two wash steps. After supernatant is discarded, carefully remove any residual liquid left at the bottom of the tube without disturbing the beads with a 10 μ l pipette tip.
- 7.10. Air dry the beads for up to 5 minutes while the tube is on the magnetic stand with the lid open.
Caution: Do not over dry the beads. This may result in lower recovery of DNA target.
- 7.11. Elute DNA target from beads by adding 20 μ l TE (0.1X) to beads. Mix well on a vortex mixer or by pipetting up and down at least 10 times. Incubate for at least 2 minutes at room temperature.
- 7.12. Quickly spin the tube and place it on a magnetic rack to separate beads from supernatant. After the solution is clear (about 2 minutes), carefully transfer 18 μ l supernatant to a new PCR tube.



Safe Stopping Point: It is safe to store the library overnight at -20°C.

8. Assess Library Quality on a Bioanalyzer High Sensitivity Chip or TapeStation High Sensitivity D1000 ScreenTape®

- 8.1 Dilute library 5-fold in nuclease-free water.
- 8.2 Run 1 μ l on a Bioanalyzer DNA High Sensitivity chip
- 8.3 Check that the electropherogram shows the expected size distribution. B Cell Receptor chain libraries will have an average size of 647 bp (Figure 4). T Cell Receptor chain libraries will have an average size of 630 bp (Figure 5). BCR + TCR chain libraries will have an average size of 632 bp (Figure 6).

Figure 4. Example of B Cell Receptor library size distribution on a Bioanalyzer.

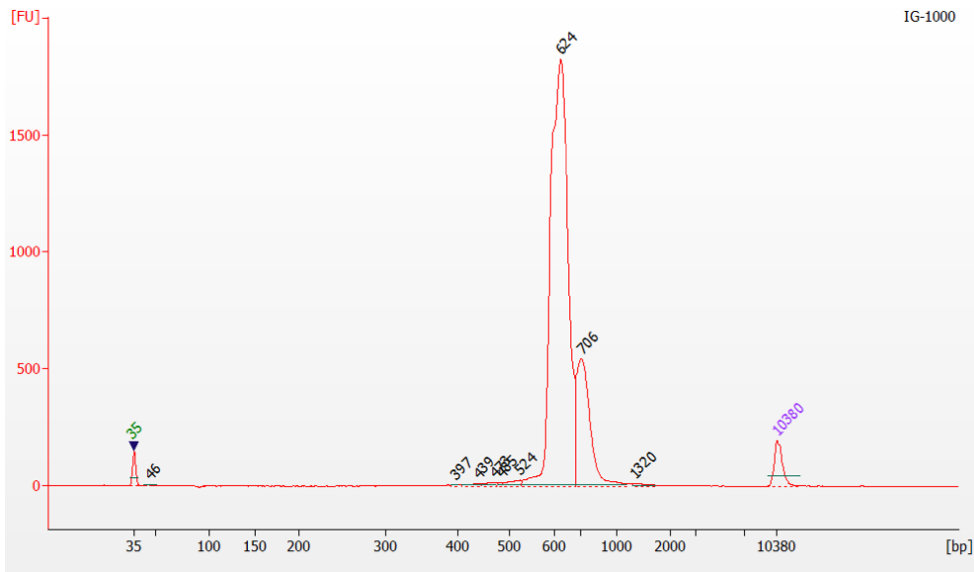


Figure 5. Example of T Cell Receptor library size distribution on a Bioanalyzer.

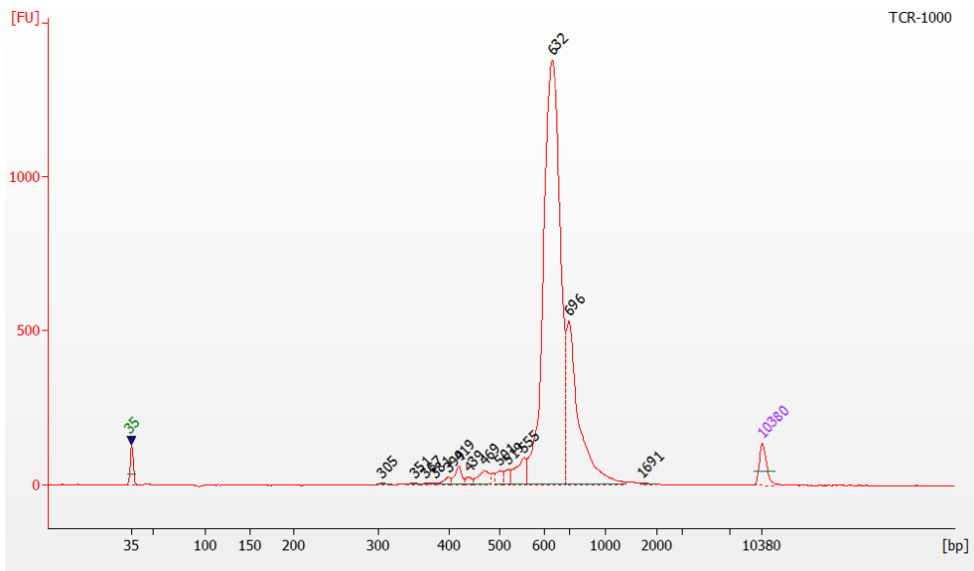
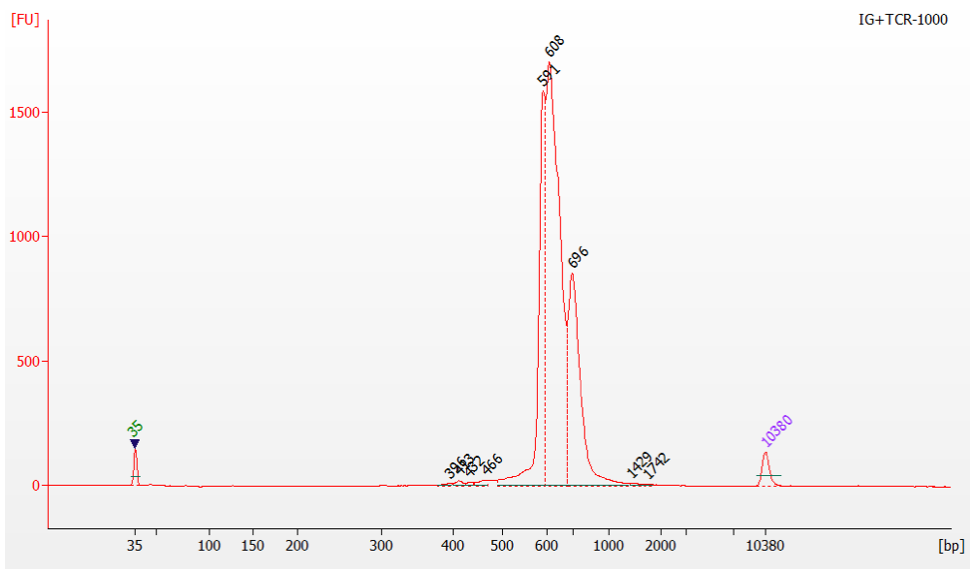


Figure 6. Example of B Cell Receptor chains and T Cell Receptor chains library size distribution on a Bioanalyzer.



9. Sequencing and Data Analysis

Sequencing should be performed on an Illumina MiSeq, using Illumina V3 sequencing chemistry. We recommend running 300 PE using a 600-cycle V3 MiSeq reagent kit with the index read between the first and second read.

For information on data processing please refer to Data Usage Guideline Page within the "Usage Guidelines & Tips" panel in the "Other Tools & Resources" tab at: www.neb.com/E6320

We have developed and optimized a pipeline for processing of data directly from FASTQ files using open-source bioinformatics tools. Details on this pipeline can be found at: <https://usegalaxy.org/u/bradlanghorst/w/presto-nebnext-immune-seq-workflow-v320>.

10. Index Pooling Guidelines

For more detailed indexing information please refer to the manual for NEBNext Multiplex Oligos for Illumina (Dual Index Primers Set 1), NEB #E7600.

For multiplexing samples, use Table 1 for some valid index combinations.

Table 1. Examples of valid index combinations.

PLEX	i7 PRIMERS	i5 PRIMERS
2	i701 and i702 i703 and i704 i705 and i706 i707 and i708 i709 and i710 i711 and i712	Any i5 primer
3	i701, i702 and i703 i703, i704 and i705 i705, i706 and i707 i707, i708 and i709 i709, i710 and i711	Any i5 primer
4	i701, i702, i703 and i704 i703, i704, i705 and i706 i705, i706, i707 and i708 i707, i708, i709 and i710 i709, i710, i711 and i712	Any i5 primer
5-12	Any valid i7 4-plex with any other i7 primers	Any i5 primer
7-12	Any 3 plex combination with and any other i7 primer (as needed)	i501, i502 and any other i5 primer (as needed) i503, i504 and any other i5 primer (as needed) i505, i506 and any other i5 primer (as needed) i507, i508 and any other i5 primer (as needed)
Greater than 12	Any 4 plex combination with any other i7 primer (as needed)	i501, i502 and any other i5 primer (as needed) i503, i504 and any other i5 primer (as needed) i505, i506 and any other i5 primer (as needed) i507, i508 and any other i5 primer (as needed)

Table 2. Index sequences of NEBNext i501 Primer – NEBNext i508 Primer:

PRODUCT	EXPECTED INDEX READ
	MiSeq
NEBNext i501 Primer	TATAGCCT
NEBNext i502 Primer	ATAGAGGC
NEBNext i503 Primer	CCTATCCT
NEBNext i504 Primer	GGCTCTGA
NEBNext i505 Primer	AGGCGAAG
NEBNext i506 Primer	TAATCTTA
NEBNext i507 Primer	CAGGACGT
NEBNext i508 Primer	GTACTGAC

Table 3. Index sequences of NEBNext i701 Primer – NEBNext i712 Primer:

PRODUCT	EXPECTED INDEX READ
NEBNext i701 Primer	ATTACTCG
NEBNext i702 Primer	TCCGGAGA
NEBNext i703 Primer	CGCTCATT
NEBNext i704 Primer	GAGATTCC
NEBNext i705 Primer	ATTCAGAA
NEBNext i706 Primer	GAATTCGT
NEBNext i707 Primer	CTGAAGCT
NEBNext i708 Primer	TAATGCGC
NEBNext i709 Primer	CGGCTATG
NEBNext i710 Primer	TCCGCGAA
NEBNext i711 Primer	TCTCGCGC
NEBNext i712 Primer	AGCGATAG

Kit Components

NEB #E6320S Table of Components

NEB #	PRODUCT	VOLUME
E6321A	NEBNext IS RT Buffer (4X)	0.120 ml
E6322A	NEBNext Cell Lysis Buffer	0.012 ml
E6323A	dNTP Solution Mix	0.108 ml
E6324A	NEBNext IS RT Primer	0.024 ml
E6326A	NEBNext IS TS Oligos	0.024 ml
E7761A	NEBNext First Strand Synthesis Enzyme Mix	0.048 ml
E6327A	Q5 Hot Start High-Fidelity DNA Polymerase	0.060 ml
E6328A	Q5 Reaction Buffer (5X)	0.600 ml
E2626A	NEBNext IS BCR Primers (Human)	0.048 ml
E2627A	NEBNext IS TCR Primers (Human)	0.048 ml
E6329A	NEBNext IS PCR2 Universal Primer	0.072 ml
E6355A	Streptavidin Magnetic Beads	0.360 ml
E6356A	NEBNext IS Bind and Wash Buffer (2X)	11.3 ml
E6357A	0.1% Tween	3.0 ml
E6358A	NEBNext Sample Purification Beads	2.8 ml
E6331A	TE (0.1X)	2.3 ml
E6332A	Nuclease-free Water	1.5 ml
E6333A	NEBNext i701 Primer	0.008 ml
E6334A	NEBNext i702 Primer	0.008 ml
E6335A	NEBNext i703 Primer	0.008 ml
E6336A	NEBNext i704 Primer	0.008 ml
E6337A	NEBNext i705 Primer	0.008 ml
E6338A	NEBNext i706 Primer	0.008 ml
E6339A	NEBNext i501 Primer	0.036 ml
E6340A	NEBNext i502 Primer	0.036 ml
E6341A	NEBNext i503 Primer	0.036 ml
E6342A	NEBNext i504 Primer	0.036 ml

NEB #E6320L Table of Components

NEB #	PRODUCT	VOLUME
E6321AA	NEBNext IS RT Buffer (4X)	0.480 ml
E6322AA	NEBNext IS Cell Lysis Buffer	0.048 ml
E6323AA	dNTP Solution Mix	0.432 ml
E6324AA	NEBNext IS RT Primer	0.096 ml
E6326AA	NEBNext IS TS Oligos	0.096 ml
E7761AA	NEBNext First Strand Synthesis Enzyme Mix	0.192 ml
E6327AA	Q5 Hot Start High-Fidelity DNA Polymerase	0.240 ml
E6328AA	Q5 Reaction Buffer (5X)	2.400 ml
E2626AA	NEBNext IS BCR Primers (Human)	0.192 ml
E6327AA	NEBNext IS TCR Primers (Human)	0.192 ml
E6329AA	NEBNext IS PCR2 Universal Primer	0.288 ml
E6355AA	Streptavidin Magnetic Beads	1.440 ml
E6356AA	NEBNext IS Bind and Wash Buffer (2X)	45.2 ml

E6357AA	0.1% Tween	11.8 ml
E6358AA	NEBNext Sample Purification Beads	11.1 ml
E6331AA	TE (0.1X)	9.2 ml
E6332AA	Nuclease-free Water	5.8 ml
E6333AA	NEBNext i701 Primer	0.032 ml
E6334AA	NEBNext i702 Primer	0.032 ml
E6335AA	NEBNext i703 Primer	0.032 ml
E6336AA	NEBNext i704 Primer	0.032 ml
E6337AA	NEBNext i705 Primer	0.032 ml
E6338AA	NEBNext i706 Primer	0.032 ml
E6347AA	NEBNext i707 Primer	0.032 ml
E6348AA	NEBNext i708 Primer	0.032 ml
E6349AA	NEBNext i709 Primer	0.032 ml
E6352AA	NEBNext i710 Primer	0.032 ml
E6353AA	NEBNext i711 Primer	0.032 ml
E6354AA	NEBNext i712 Primer	0.032 ml
E6339AA	NEBNext i501 Primer	0.144 ml
E6340AA	NEBNext i502 Primer	0.144 ml
E6341AA	NEBNext i503 Primer	0.144 ml
E6342AA	NEBNext i504 Primer	0.144 ml
E6343AA	NEBNext i505 Primer	0.144 ml
E6344AA	NEBNext i506 Primer	0.144 ml
E6345AA	NEBNext i507 Primer	0.144 ml
E6346AA	NEBNext i508 Primer	0.144 ml

CheckList

1. NEBNext Immune Sequencing Reverse Transcription and cDNA Synthesis

- 1.1. Add RT Reagents to up to 9 μ l total RNA Sample to make 20.5 μ l total reaction volume:
 - 5 μ l RT Buffer (4X)
 - 1 μ l RT Primer
 - 1 μ l TS Oligos
 - 2 μ l dNTP Solution Mix
 - 2 μ l First Strand Synthesis Enzyme Mix
 - 0.5 μ l Cell Lysis Buffer
 - Add ___ μ l Nuclease-free water
- 1.2. Pipette mix 10 times with pipette set to 15 μ l, quick spin
- 1.3. Thermal cycler (Heated lid \geq 80°C; 40 minutes 42°C, 10 minutes 70°C, Hold at 4°C)

2. Purify cDNA

- 2.1. Prepare NEBNext IS Bind and Wash Buffer (1X):
 - 50 μ l NEBNext IS Bind and Wash Buffer (2X)
 - 50 μ l nuclease-free water
- 2.2. Vortex Streptavidin Magnetic Beads and add 15 μ l to a new 1.5 ml tube
- 2.3. Place tube on magnet, wait 2 min and remove supernatant
- 2.4. 200 μ l undiluted NEBNext IS Bind and Wash Buffer (2X), vortex and quick spin
- 2.5. Place tube on magnet, wait until solution is clear and remove supernatant
- 2.6. Repeat Step 2.4–2.5 once
- 2.7. Add 20 μ l undiluted NEBNext IS Bind and Wash Buffer (2X) to beads, vortex and quick spin
- 2.8. Add 20 μ l beads to cDNA product, vortex and quick spin
- 2.9. Rotate at room temperature for 15 min
- 2.10. Quick spin, place on magnet, wait until solution is clear and remove supernatant
- 2.11. Add 100 μ l diluted NEBNext IS Bind and Wash Buffer (1X)
- 2.12. Quick spin, place on magnet, wait until solution is clear and remove supernatant
- 2.13. Add 100 μ l 0.1% Tween and vortex
- 2.14. Quick spin, place on magnet, wait until solution is clear and remove supernatant
- 2.15. Add 23 μ l 0.1% Tween and vortex and quick spin
- 2.16. Thermal cycler (Heated lid \geq 105°C; 3 minutes 95°C, Hold at 25°C)
- 2.17. Vortex, quick spin and place on magnet
- 2.18. Wait until solution is clear and transfer 21 μ l to a new tube

3. PCR1

- 3.1. Add PCR1 reagents to 21 μ l cDNA sample. (Choose between steps 3.1A, 3.1B and 3.1C):
 - 3.1A. PCR1 for Enriching B Cell Receptor Chains
 - 10 μ l Q5 Reaction Buffer (5X)
 - 1 μ l dNTP Solution Mix
 - 1 μ l Q5 Hot Start High-Fidelity DNA Polymerase
 - 2 μ l NEBNext IS BCR Primers (Human)
 - 1 μ l NEBNext i7 Index Primer
 - 14 μ l Nuclease-free Water

- 3.1B. PCR1 for Enriching T Cell Receptor Chains
- 10 µl Q5 Reaction Buffer (5X)
 - 1 µl dNTP Solution Mix
 - 1 µl Q5 Hot Start High-Fidelity DNA Polymerase
 - 2 µl NEBNext IS TCR Primers (Human)
 - 1 µl NEBNext i7 Index Primer
 - 14 µl Nuclease-free Water
- 3.1C. PCR1 for Enriching Both B Cell Receptor Chains and T Cell Receptor Chains
- 10 µl Q5 Reaction Buffer (5X)
 - 1 µl dNTP Mix
 - 1 µl Q5 Hot Start High-Fidelity DNA Polymerase
 - 2 µl NEBNext IS BCR Primers (Human)
 - 2 µl NEBNext IS TCR Primers (Human)
 - 1 µl NEBNext i7 Index Primer
 - 12 µl Nuclease-free Water
- 3.2. Pipette mix 10 times with pipette set to 40 µl, quick spin
- 3.3. Thermal cycler (Heated lid $\geq 105^{\circ}\text{C}$):
- 60 seconds at 98°C ,
 - 12 cycles of (10 seconds at 98°C , 30 seconds at 64°C , 30 seconds at 72°C)
 - 60 seconds at 72°C
 - Hold at 4°C

4. Cleanup of PCR1 Amplification

- 4.1. Vortex NEBNext Sample Purification Beads
- 4.2. Add 50 µl beads to sample and mix by pipetting 10 times
- 4.3. Incubate 5 min
- 4.4. Place tubes on magnet, wait 5 min then remove the supernatant (keep the beads)
- 4.5. On magnet add 200 µl 80% ethanol, wait 30 seconds and remove
- 4.6. Repeat Step 4.5. once
- 4.7. Wait 5 min with tubes on magnet and lid open
- 4.8. Add 25 µl TE (0.1X) and mix by pipetting 10 times
- 4.9. Place tubes on magnet, wait 2 min then transfer 23 µl to new tubes

5. qPCR for PCR2 Cycle Optimization

- 5.1. Add qPCR reagents to 5 µl of the sample:
 - 5 µl Q5 Reaction Buffer (5X)
 - 0.5 µl dNTP Solution Mix
 - 0.5 µl Q5 Hot Start High-Fidelity DNA Polymerase
 - 1 µl NEBNext IS PCR2 Universal Primer
 - 1 µl NEBNext i5 Primer
 - 0.25 µl SYBR Green (20X)
 - 11.75 µl Nuclease-free Water
- 5.2. Pipette mix 10 times with pipette set to 20 µl, quick spin

5.3. qPCR program:
60 seconds at 98°C
30 cycles of (10 seconds at 98°C; 30 seconds at 72°C, fluorescence on)
Hold at 4°C

5.4. Determine cycle number at 2/3 of peak fluorescence

6. PCR2 for Final Library

6.1. Add PCR2 reagents to the remaining 10 µl of PCR sample:

10 µl Q5 Reaction Buffer (5X)

1 µl dNTP Solution Mix

1 µl Q5 Hot Start High-Fidelity DNA Polymerase

2 µl NEBNext IS PCR2 Universal Primer

2 µl NEBNext i5 Primer

24 µl Nuclease-free Water

6.2. Pipette mix 10 times with pipette set to 40 µl, quick spin

6.3. Thermal cycler (Heated lid \geq 105°C):

60 seconds at 98°C

__ cycles of (10 seconds at 98°C; 30 seconds at 72°C)

Hold at 4°C

7. Cleanup of PCR2 Amplification

7.1. Vortex NEBNext Sample Purification Beads

7.2. Add 50 µl TE (0.1X) and 50 µl beads to sample and mix by pipetting 10 times

7.3. Incubate 5 min

7.4. Place tubes on magnet, wait 5 min then then transfer the supernatant to a new tube (keep the supernatant)

7.5. Add 15 µl NEBNext Sample Purification Beads to sample and mix by pipetting 10 times

7.6. Incubate 5 min

7.7. Place tubes on magnet, wait 5 min then remove the supernatant (keep the beads)

7.8. On magnet add 200 µl 80% ethanol, wait 30 seconds and remove

7.9. Repeat Step 7.8. once

7.10. Wait 5 min with tubes on magnet and lid open

7.11. Add 20 µl TE (0.1X) and mix by pipetting 10 times

7.12. Place tubes on magnet, wait 2 min then transfer 18 µl to new tubes

8. Library QC on Bioanalyzer

8.1. Dilute library 5-fold in nuclease-free water

8.2. Run 1µl on a Bioanalyzer High Sensitivity DNA chip

8.3. Assess library quality

Revision History

REVISION #	DESCRIPTION	DATE
1.0		N/A
2.0	Update protocol text.	12/21

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

This product is covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information about commercial rights, please email us at busdev@neb.com. While NEB develops and validates its products for various applications, the use of this product may require the buyer to obtain additional third party intellectual property rights for certain applications.

This product is licensed for research and commercial use from Bio-Rad Laboratories, Inc., under U.S. Pat. Nos. 6,627,424, 7,541,170, 7,670,808, 7,666,645, and corresponding patents in other countries. No rights are granted for use of the product for Digital PCR or real-time PCR applications, with the exception of quantification in Next Generation Sequencing workflows.

This product is licensed for preclinical and clinical research uses, and commercial applications related to the foregoing, but excluding: (a) immune sequencing services (i.e., performance of immune sequencing services on a for-fee basis or otherwise in exchange for consideration); (b) immune sequencing partnerships; (c) immune sequencing collaborations; (d) research, development or commercialization of FDA [or other regulatory authority outside the U.S. having similar regulatory authority as the FDA for such other jurisdiction(s)]-approved diagnostic products; and (e) any *in vitro* diagnostic products.

AMPURE® are registered trademarks of Beckman Coulter, Inc.

B CORPORATION™ is a trademark of B Lab, Inc.

ILLUMINA® HISEQ®, MISEQ®, NEXTSEQ®, NOVASEQ® and MINISEQ® are registered trademarks of Illumina, Inc.

BIOANALYZER®, SCREENTAPE® and TAPESTATION® are registered trademarks of Agilent Technologies

SYBR® is a registered trademark of Molecular Probes, Inc.

LOBIND® is a registered trademark of Eppendorf AG.

© Copyright 2021, New England Biolabs, Inc.; all rights reserved



www.neb.com



be INSPIRED
drive DISCOVERY
stay GENUINE