

NEBNext® ARTIC SARS-CoV-2 Companion Kit (Oxford Nanopore Technologies®)

NEB #E7660S/L

24/96 reactions

Version 8.0_5/26

Table of Contents

Overview	2
Chapter 1	
NEBNext ARTIC Express Protocol without PCR Bead Cleanup (Two clean-up steps)	4
Chapter 2	
NEBNext ARTIC Standard Protocol with PCR Bead Cleanup (Three clean-up steps)	10
Chapter 3	
NEBNext VarSkip Short v2 Express Protocol without PCR Bead Cleanup (Two clean-up steps)	17
Kit Components	23
Revision History	24

The Kit Includes

The volumes provided are sufficient for preparation of up to 24 barcoding reactions (NEB #E7660S, minimum 6 barcoding samples per run for total 4 runs) and 96 barcoding reactions (NEB #E7660L, minimum 24 barcoding samples per run for total 4 runs). Colored bullets represent the color of the cap of the tube containing the reagent. If one plans to follow a different protocol, additional reagents can be purchased separately)

Package 1: Store at –20°C. (NEB #E7660-1, 24 reactions, NEB #E7660-2, 96 reactions)

- (lilac) LunaScript® RT SuperMix
- (lilac) Q5® Hot Start High-Fidelity 2X Master Mix
- (green) NEBNext Ultra II End Prep Enzyme Mix
- (green) NEBNext Ultra II End Prep Reaction Buffer
- (red) Blunt/TA Ligase Master Mix
- (red) NEBNext Quick T4 DNA Ligase
- (red) NEBNext Quick Ligation Reaction Buffer
- (white) Nuclease-free water

Package 2: Store at –20°C. (NEB #E7871S, 24 reactions and NEB #E7871L, 96 reactions)

- (lilac) NEBNext ARTIC SARS-CoV-2 Primer Mix 1
- (lilac) NEBNext ARTIC SARS-CoV-2 Primer Mix 2
- (lilac) NEBNext ARTIC Human Control Primer Pairs 1
- (lilac) NEBNext ARTIC Human Control Primer Pairs 2
- (orange) NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1
- (orange) NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2

Package 3: Store at room temperature. Do not freeze. (NEB #E7666S, 24 reactions and NEB #E7666L, 96 reactions)

NEBNext Sample Purification Beads

Required Materials Not Included

- 80% Ethanol (freshly prepared)
- DNA LoBind Tubes (Eppendorf® #022431021)
- Oxford Nanopore Technologies Native Barcoding Expansion Kits 1-12 (EXP-NBD104), 13-24 (EXP-NBD114) or 1-96 (EXP-NBD196)
- Oxford Nanopore Technologies Ligation sequencing DNA V14 (SQK-LSK114) or Oxford Nanopore Technologies Adapter Mix II Expansion (EXP-AMII001), Flow Cell Priming kit (EXP-FLP002) and Sequencing Auxiliary Vials (EXP-AUX001)
- Oxford Nanopore Technologies SFB Expansion Kit (EXP-SFB001)
- Qubit® dsDNA HS Assay Kit (Thermo Fisher Scientific, Inc.® Q32851)
- Magnetic rack/stand (NEB #S1515, Alpaqua®, cat. #A001322 or equivalent)
- Thermal cycler
- Vortex Mixer
- Microcentrifuge
- Agilent® Bioanalyzer® or similar fragment analyzer and associated consumables (#4150 or #4200 TapeStation System)
- DNase RNase free PCR strip tubes (USA Scientific 1402-1708)
- 1.5 ml tube magnet stand (NEB #S1506)

Protocol Descriptions

Chapter 1: NEBNext ARTIC Express Protocol without PCR Bead Cleanup (Two clean-up steps): This protocol utilizes NEBNext ARTIC SARS-CoV-2 Primer Mixes for targeting SARS-CoV-2, these are balanced ARTICv3 primers, This protocol does not include a cleanup and normalization step for each sample after cDNA synthesis. Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. Skipping these steps reduces hands on time but may require a longer sequencing run to obtain sufficient coverage for each sample.

Chapter 2: NEBNext ARTIC Standard Protocol with PCR Bead Cleanup (Three clean-up steps): This protocol includes a cleanup and normalization step for each sample after cDNA synthesis. Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. These pools will likely achieve sufficient and equal coverage in less run time, but they take more hands-on time.

Protocols.io also provides an interactive version of this protocol where you can discover and share optimizations with the research community. Please look for E7660 on <https://www.protocols.io/>

Chapter 3: NEBNext VarSkip Short v2 Express Protocol without PCR Bead Cleanup (Two clean-up steps): This protocol follows an alternate variant-tolerant approach for targeting SARS-CoV-2 by utilizing NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mixes. The NEBNext VarSkip Short v2 SARS-CoV-2 Primer mixes cannot be added to the same cDNA amplification reaction as the NEBNext ARTIC SARS-CoV-2 Primer Mixes. This protocol does not include a cleanup and normalization step for each sample after cDNA synthesis. Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. Skipping these steps reduces hands on time but may require a longer sequencing run to obtain sufficient coverage for each sample.

Chapter 1

NEBNext ARTIC Express Protocol without PCR Bead Cleanup (Two clean-up steps)

This protocol does not include a cleanup and normalization step for each sample after cDNA synthesis. Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. Skipping these steps reduces hands on time, but may require a longer sequencing run to obtain sufficient coverage for each sample. To obtain more even sample to sample coverage, we recommend normalizing the RNA samples prior to starting Express Protocol without PCR Bead Cleanup (this Chapter).

Symbols



This is a point where you can safely stop the protocol.



This caution sign signifies a step in the protocol that has two paths leading to the same end point but is dependent on a user variable, like the amount of input DNA.



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

Note 1: The amount of RNA required for detection depends on the abundance of the RNA of interest. In general, we recommend, using ≥ 10 copies of the SARS-CoV-2 viral genome as input. The presence of carry-over products can interfere with sequencing accuracy, particularly for low copy targets. Therefore, it is important to carry out the appropriate no template control (NTC) reactions to demonstrate that positive reactions are meaningful. We recommend that the NTC reactions are prepared prior to non-NTC samples following the guidelines outlined below.

Guidelines for Sample Setup: We recommend setting up the cDNA synthesis and cDNA PCR setup reactions in a room (and ideally in a hood) separate from cDNA amplification and library construction area to minimize cross-contamination of future reactions.

Note 2: If sample Ct is between 12-15, then it is recommended per the ARTIC network [nCoV 2019 sequencing protocol v3 LoCost](#) to dilute the sample 100-fold in water, if between 15-18 then dilute 10-fold in water. This will reduce the likelihood of PCR-inhibition.

1.1. cDNA Synthesis

1.1.1. Gently mix 10 times by pipetting and spin down the LunaScript RT SuperMix reagents (contains primers). Prepare the cDNA synthesis reaction as described below:

COMPONENT	VOLUME
RNA Sample*	8 μ l
• (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

*Up to 0.5 μ g total RNA can be used in a 10 μ l reaction.

1.1.2. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

1.1.3. For no template controls, mix the following components

COMPONENT	VOLUME
○ (white) Nuclease-free Water	8 μ l
• (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

1.1.4. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

1.1.5. Incubate reactions in a thermal cycler with lid temperature at 105°C with the following steps:

CYCLE STEP	TEMP	TIME	CYCLE
Primer Annealing	25°C	2 minutes	1
cDNA Synthesis	55°C	20 minutes	
Heat Inactivation	95°C	1 minute	

*Set heated lid to 105°C



Samples can be stored at –20°C if they are not used immediately.

1.2. Targeted cDNA Amplification

Note: 4.5 µl cDNA input is recommended. If using less than 4.5 µl of cDNA, add nuclease-free water to a final volume of 4.5 µl.

Use of the NEBNext ARTIC Human Control Primer Pairs 1 and 2 are optional. If used, the appropriate NEBNext ARTIC Human Control Primer Pairs and NEBNext ARTIC SARS-CoV-2 Primer Mix should be combined prior to use. More specifically, NEBNext ARTIC Human Control Primer Pairs 1 should be combined with NEBNext ARTIC SARS-CoV-2 Primer Mix 1 and NEBNext ARTIC Human Control Primer Pairs 2 with NEBNext ARTIC SARS-CoV-2 Primer Mix 2. Mixing directions are listed below.

- 1.2.1. Gently mix Q5 Hot Start High Fidelity 2X Master Mix 10 times by pipetting and spin down reagents. Prepare the split pool amplification reactions as described below.

For Pool Set A:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 1 with 42 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 1, vortex and spin down reagents. If using a 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Mix 1 with 168 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 1, vortex and spin down reagents. Use 1.75 µl of the combined mix for each Pool Set A reaction.

COMPONENT	VOLUME
cDNA (Step 1.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext ARTIC SARS-CoV-2 Primer Mix 1*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 1, add 1.75 µl of the combined NEBNext ARTIC Human Control Primer Pairs 1 and NEBNext ARTIC SARS-CoV-2 Primer Mix 1.

- 1.2.2. Flick the tube or pipet up and down to mix followed by a quick spin.

1.2.3. For Pool Set B:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 2 with 42 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 2, vortex and spin down reagents. If using 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Pairs 2 with 168 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 2. Use 1.75 µl of the combined mix for each Pool Set B reaction.

COMPONENT	VOLUME
cDNA (Step 1.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext ARTIC SARS-CoV-2 Primer Mix 2*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 2, add 1.75 µl of the combined NEBNext ARTIC Human Control Primer Pairs 2 and NEBNext ARTIC SARS-CoV2 Primer Mix 2.

Note: We recommend proceeding with the cDNA amplification and library construction in a different area or room.

- 1.2.4. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

- 1.2.5. Incubate reactions in a thermal cycler** with the following steps:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	30 second	1
Denature	95°C	15 seconds	35
Annealing and Extension	63°C*	5 minutes	
Hold	4°C	∞	1

* It is very important to set up the annealing and extension temperature to 63°C

** Set heated lid to 105°C.



Samples can be stored at –20°C if they are not used immediately.

1.3. PCR Reaction Pooling

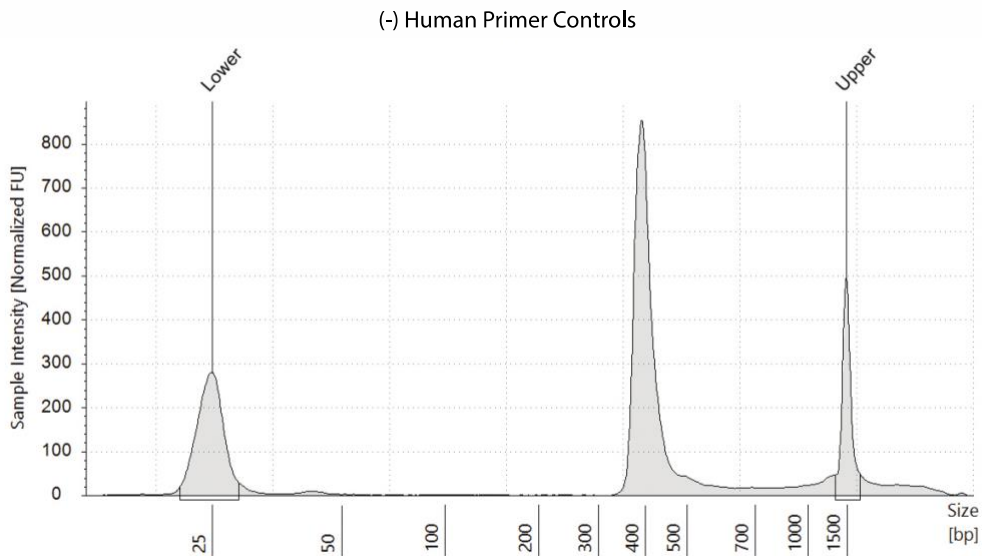
Please note, there is also a protocol that includes a cleanup and normalization step at this point (chapter 2 in this manual). Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. These pools will likely achieve sufficient coverage in less run time

- 1.3.1. For each sample, combine pool A and pool B PCR Reactions.

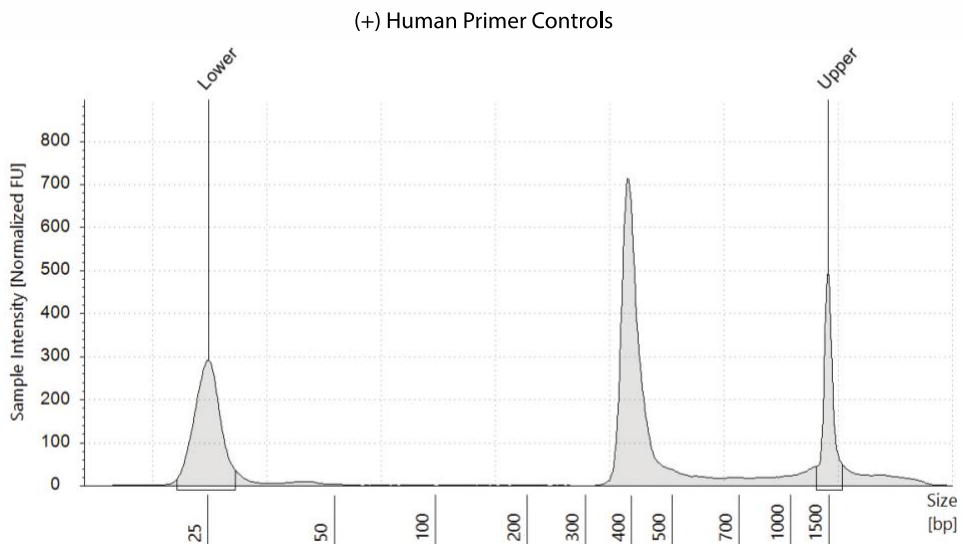


Figure 1.3.1: Example of cDNA amplicons generated from 1000 genome copies of SARS CoV-2 in the absence (A) and presence (B) of human primer controls.

A



B



1.4. NEBNext End Prep

- 1.4.1. Add the following components to a PCR tube (End Prep Reaction and Buffer can be pre-mixed and master mix is stable on ice for 4 hours):

COMPONENT	VOLUME
Targeted cDNA Amplicons (1.3.1)	1 µl
○ (white) Nuclease-free water	11.5 µl
● (green) NEBNext Ultra II End Prep Reaction Buffer	1.75 µl
● (green) NEBNext Ultra II End Prep Enzyme Mix	0.75 µl
Total Volume	15 µl

- 1.4.2. Flick the tube or pipet up and down 10 times to mix the solution. 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 performance.

- 1.4.3. Place in a thermal cycler, with the heated lid set to $\geq 75^{\circ}\text{C}$, and run the following program:

10 minutes @ 20°C

10 minutes @ 65°C

Hold at 4°C



If necessary, samples can be stored at -20°C for a few days; however, a slight loss in yield (~20%) may be observed. We recommend continuing with barcode ligation before stopping.

1.5. Barcode Ligation

- 1.5.1. Add the following components directly to a sterile nuclease-free PCR tube:

COMPONENT	VOLUME
○ (white) Nuclease-free water	6 µl
End-prepped DNA (1.4.3.)	1.5 µl
Native Barcode*	2.5 µl
● (red) Blunt/TA Ligase Master Mix**	10 µl
Total Volume	20 µl

* Native Barcodes are provided in Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104) and 13-24 (EXP-NBD114) or 1-96 (EXP-NBD196)

** Mix the Blunt/TA Ligase Master Mix by pipetting up and down several times prior to adding to the reaction

- 1.5.2. Flick the tube or pipet up and down 10 times to mix solution. Perform a quick spin to collect all liquid from the sides of the tube. **(Caution: The Blunt/TA Ligase Master Mix is very viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**

- 1.5.3. Incubate at 25°C for 20 minutes.
Incubate at 65°C for 10 minutes.
Place on ice for 1 min.

- 1.5.4. Pool all barcoded samples into one 1.5 ml DNA LoBind Tube.

1.6. Cleanup of Barcoded DNA

Note: Use the pooled barcoded DNA samples (from Step 1.5.4.), up to 480 µl for bead cleanup. Remaining pooled DNA can be stored at -20°C

- 1.6.1. Vortex NEBNext Sample Purification Beads to resuspend.

- 1.6.2. Add 0.4X resuspended beads to pooled barcoded samples (1.5.4.) (for example, if you are pooling 24 barcoded DNA samples [which amounts to 480 µl total], add 192 µl of resuspended sample purification beads to the 480 µl of pooled sample). Mix well by flicking the tube or pipetting up and down 10 times. Perform a quick spin for 1 second to collect all liquid from the sides of the tube.
- 1.6.3. Incubate samples on bench top for 10 minutes at room temperature.
- 1.6.4. Place the tube on a 1.5 ml magnetic stand (such as NEB #S1506) to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.
- 1.6.5. After 3 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard beads**).
- 1.6.6. Wash the beads by adding 250 µl of Short Fragment buffer (SFB). Flick the tube or pipet up and down 10 times to mix to resuspend pellet. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 1.6.7. Place the tube on an appropriate magnetic stand for 3 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.
- 1.6.8. Repeat step 1.6.6. and 1.6.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube, place back on the magnetic stand and remove traces of SFB with a p10 pipette tip
- 1.6.9. Add 500 µl of 80% freshly prepared ethanol to the tube while on the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets. Repeat the wash once to make it a total of two washes.
- 1.6.10. A quick spin and place the sample tube on the magnetic stand, remove any residual ethanol.
- 1.6.11. Air dry the beads for 30 seconds 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. Elute the samples when the beads are still dark brown and glossy looking, but when all visible liquid has evaporated. When the beads turn lighter brown and start to crack, they are too dry.**
- 1.6.12. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 33 µl of Nuclease-free water.
- 1.6.13. Resuspend the pellet by flicking the tube or pipetting up and down 10 times to mix. Incubate for at least 2 minutes at room temperature. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube before placing back on the magnetic stand.
- 1.6.14. Place the tube on the magnetic stand. After 2 minutes (or when the solution is clear), transfer 32 µl to a new 1.5 ml Eppendorf DNA LoBind Tube or PCR tube.
- 1.6.15. Assess the concentration of the purified barcoded DNA sample. We recommend using a Qubit fluorometer for concentration assessment. Nanodrop is NOT recommended since it may overestimate the DNA concentration. Use 1 µl of sample for the Qubit fluorometer.



Samples can be stored at –20°C if they are not used immediately.

1.7. Adapter Ligation

- 1.7.1. Use the Qubit readings from Step 1.6.15. to dilute up to 60 ng of the Native barcoded DNA pool with nuclease-free water to a final volume of 30 µl (2 ng/µl). Add the following components into a 1.5 ml Eppendorf DNA LoBind Tube or nuclease-free PCR tube:

COMPONENT	VOLUME
Native barcoded and purified DNA (1.6.14.; up to 60 ng)	30 µl
● (red) NEBNext Quick Ligation Reaction Buffer*	10 µl
Adapter Mix II (AMII)**	5 µl
● (red) NEBNext Quick T4 Ligase	5 µl
Total Volume	50 µl

* Mix the NEBNext Quick Ligation Reaction Buffer by pipetting up and down several times prior to adding to the reaction.

** Adapter Mix II is provided by Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104), 13-24 (EXP-NBD114) and 1-96 (EXP-NBD-196) kits.

- 1.7.2. Flick the tube to mix solution. Perform a quick spin for 1 second to collect all liquid from the sides of the tube. **(Caution: The NEBNext Quick Ligation Buffer is viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**
- 1.7.3. Incubate at 25°C or room temperature for 20 minutes.
- 1.7.4. Proceed to Cleanup of Adapter-ligated DNA in Section 1.8.

1.8. Cleanup of Adapter Ligated DNA

Note: The volumes of SPRIselect or NEBNext Sample Purification Beads provided here are for use with the sample contained in the exact buffer at this step. AMPure XP beads can be used as well. If using AMPure XP beads, allow the beads to warm to room temperature for at least 30 minutes before use. These volumes may not work properly for a cleanup at a different step in the workflow.

- 1.8.1. Vortex NEBNext Sample Purification Beads to resuspend.
- 1.8.2. Add 50 µl (1X) resuspended beads to the ligation mix. Mix well by flicking the tube followed by a quick spin for 1 second.
- 1.8.3. Incubate samples for 10 minutes at room temperature.
- 1.8.4. Place the tube on an appropriate magnetic stand to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.
- 1.8.5. After 3 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets **(Caution: do not discard the beads)**.
- 1.8.6. Wash the beads by adding 250 µl of Short Fragment Buffer (SFB). Flick the tube to resuspend pellet. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand. Place the tube on an appropriate magnetic stand.
- 1.8.7. Wait for 3 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.
- 1.8.8. Repeat steps 1.8.6. and 1.8.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube/plate, place back on the magnet and remove traces of SFB with a p10 pipette tip.
- 1.8.9. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 15 µl of Elution Buffer (EB) provided in SQK-LSK114 kit from Oxford Nanopore.
- 1.8.10. Resuspend the pellet well in EB buffer by flicking. Incubate for 10 minutes at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 1.8.11. Place the tube/plate on the magnetic stand. After 3 minutes (or when the solution is clear), transfer 15 µl to a new DNA LoBind tube.
- 1.8.12. Use Qubit to quantify 1 µl DNA sample. Follow Oxford Nanopore Protocol SQK-LSK114 to prepare MinION® flow cell and DNA library sequencing mix using up to 20 ng adapter-ligated DNA sample (Step 1.8.11).

Note: After normalizing the DNA to 20 ng, if the volume is less than 12 µl, then top off the sample volume to 12 µl with EB.

Chapter 2

NEBNext ARTIC Standard Protocol with PCR Bead Cleanup (Three clean-up steps)

This protocol includes a cleanup and normalization step for each sample after cDNA synthesis. Performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. These pools will likely achieve sufficient and equal coverage in less run time, but they take more hands-on time.

Symbols



This is a point where you can safely stop the protocol.



This caution sign signifies a step in the protocol that has two paths leading to the same end point but is dependent on a user variable, like the amount of input DNA.



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

Note 1: The amount of RNA required for detection depends on the abundance of the RNA of interest. In general, we recommend, using ≥ 10 copies of the SARS-CoV-2 viral genome as input. The presence of carry-over products can interfere with sequencing accuracy, particularly for low copy targets. Therefore, it is important to carry out the appropriate no template control (NTC) reactions to demonstrate that positive reactions are meaningful. We recommend that the NTC reactions are prepared prior to non-NTC samples following the guidelines outlined below.

Guidelines for Sample Setup: We recommend setting up the cDNA synthesis and cDNA PCR setup reactions in a room (and ideally in a hood) separate from cDNA amplification and library construction area to minimize cross-contamination of future reactions.

Note 2: If sample Ct is between 12-15, then it is recommended per the ARTIC network [nCoV 2019 sequencing protocol v3 LoCost](#) to dilute the sample 100-fold in water, if between 15-18 then dilute 10-fold in water. This will reduce the likelihood of PCR-inhibition.

2.1. cDNA Synthesis

2.1.1. Gently mix 10 times by pipetting and spin down the LunaScript RT SuperMix reagents (contains primers). Prepare the cDNA synthesis reaction as described below:

COMPONENT	VOLUME
RNA Sample*	8 μ l
• (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

*Up to 0.5 μ g total RNA can be used in a 10 μ l reaction.

2.1.2. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

2.1.3. For no template controls, mix the following components

COMPONENT	VOLUME
○ (white) Nuclease-free Water	8 μ l
• (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

2.1.4. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

2.1.5. Incubate reactions in a thermal cycler with lid temperature at 105°C with the following steps:

CYCLE STEP	TEMP	TIME	CYCLE
Primer Annealing	25°C	2 minutes	1
cDNA Synthesis	55°C	20 minutes	
Heat Inactivation	95°C	1 minute	



Samples can be stored at –20°C if they are not used immediately.

2.2. Targeted cDNA Amplification

Note: 4.5 µl cDNA input is recommended. If using less than 4.5 µl of cDNA, add nuclease-free water to a final volume of 4.5 µl.

Use of the NEBNext ARTIC Human Control Primer Pairs 1 and 2 are optional. If used, the appropriate NEBNext ARTIC Human Control Primer Pairs and NEBNext ARTIC SARS-CoV-2 Primer Mix should be combined prior to use. More specifically, NEBNext ARTIC Human Control Primer Pairs 1 should be combined with NEBNext ARTIC SARS-CoV-2 Primer Mix 1 and NEBNext ARTIC Human Control Primer Pairs 2 with NEBNext ARTIC SARS-CoV-2 Primer Mix 2. Mixing directions are listed below.

- 2.2.1. Gently mix Q5 Hot Start High Fidelity 2X Master Mix 10 times by pipetting and spin down reagents. Prepare the split pool amplification reactions as described below.

For Pool Set A:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 1 with 42 µl of ARTIC SARS-CoV-2 Primer Mix 1, vortex and spin down reagents. If using a 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Mix 1 with 168 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 1, vortex and spin down reagents. Use 1.75 µl of the combined mix for each Pool Set A reaction.

COMPONENT	VOLUME
cDNA (Step 2.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext ARTIC SARS-CoV-2 Primer Mix 1*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 1, add 1.75 µl of the combined NEBNext ARTIC Human Control Primer Pairs 1 and NEBNext ARTIC SARS-CoV-2 Primer Mix 1.

- 2.2.2. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

2.2.3. For Pool Set B:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 2 with 42 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 2, vortex and spin down reagents. If using 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Pairs 2 with 168 µl of NEBNext ARTIC SARS-CoV-2 Primer Mix 2. Use 1.75 µl of the combined mix for each Pool Set B reaction.

COMPONENT	VOLUME
cDNA (Step 2.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext ARTIC SARS-CoV-2 Primer Mix 2*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 2, add 1.75 µl of the combined NEBNext ARTIC Human Control Primer Pairs 2 and NEBNext ARTIC SARS-CoV-2 Primer Mix 2.

Note: We recommend proceeding with the cDNA amplification and library construction in a different area or room.

- 2.2.4. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

- 2.2.5. Incubate reactions in a thermal cycler* with the following steps:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	30 second	1
Denature	95°C	15 seconds	35
Annealing and Extension	63°C**	5 minutes	
Hold	4°C	∞	1

* Set heated lid to 105°C.

** It is very important to set up the annealing and extension temperature to 63°C



Samples can be stored at –20°C if they are not used immediately.

2.3. Cleanup of cDNA Amplicons

If you prefer to omit the cleanup follow chapter 1 in this manual. **Otherwise continue to Cleanup of cDNA Amplicons (Step 2.3.1).**



Note: SPRIselect or AMPure® XP Beads can be used as well. If using AMPure XP Beads, allow the beads to warm to room temperature for at least 30 minutes before use. These bead volumes may not work properly for a cleanup at a different step in the workflow. For cleanups of samples contained in different buffer conditions, the volumes may need to be experimentally determined.

- 2.3.1. For each sample, combine pool A and pool B PCR Reactions.
- 2.3.2. Vortex SPRIselect or NEBNext Sample Purification Beads to resuspend.
- 2.3.3. Add 20 µl (0.8X) resuspended beads to the combined PCR reaction. Mix well by flicking the tube or pipetting up and down 10 times to mix and a very short 2-3 seconds quick centrifugation. Be sure to stop the centrifugation before the beads start to settle out.
- 2.3.4. Incubate samples at room temperature for 10 minutes.
- 2.3.5. Place the tubes on an appropriate magnetic stand to separate the beads from the supernatant. If necessary, quickly spin the sample 1 second to collect the liquid from the sides of the tube before placing on the magnetic stand.
- 2.3.6. After 2 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard the beads**).
- 2.3.7. Add 500 µl of 80% freshly prepared ethanol to the tube while in the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets.
- 2.3.8. Repeat step 2.3.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube for 1 second, place back on the magnetic stand and remove traces of ethanol with a p10 pipette tip.
- 2.3.9. Air dry the beads for 30 seconds 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. Elute the samples when the beads are still dark brown and glossy looking. When the beads turn lighter brown and start to crack, they are too dry.

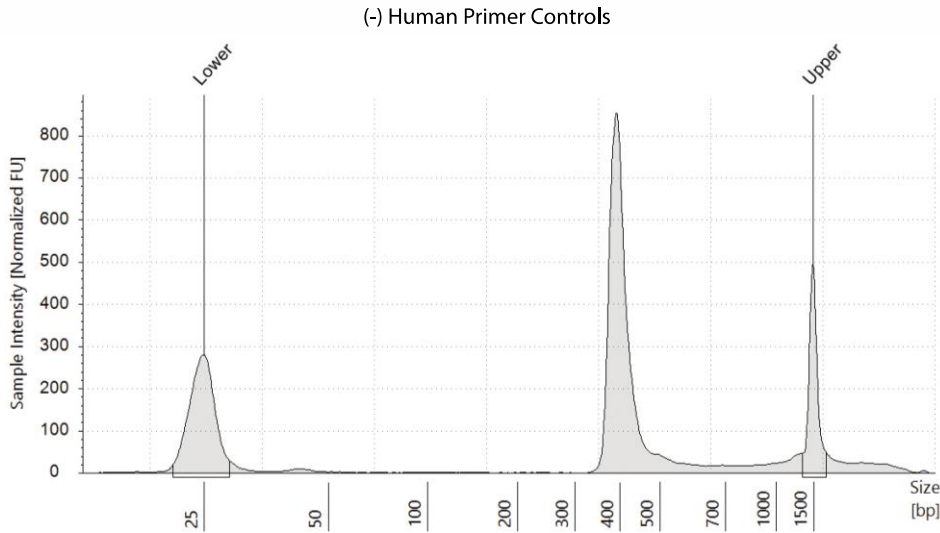
- 2.3.10. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 21 µl Nuclease-free water.
- 2.3.11. Mix well by flicking the tube or pipetting up and down 10 times to mix and followed by a very short centrifugation. Incubate for 10 minutes at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 2.3.12. Place the tube on the magnetic stand. After 2 minutes (or when the solution is clear), transfer 20 µl to clean PCR tubes.
- 2.3.13. Assess the concentration of the purified samples. We recommend using a Qubit fluorometer for concentration assessment. Use 1 µl of sample for the Qubit fluorometer.



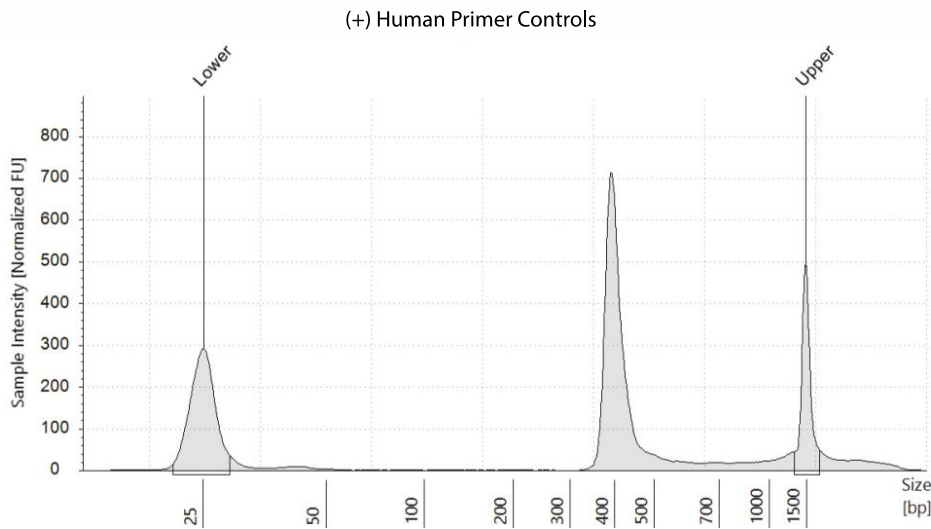
Samples can be stored at –20°C if they are not used immediately.

Figure 2.3.13: Example of cDNA amplicons generated from 1000 genome copies of SARS CoV-2 in the absence (A) and presence (B) of human primer controls.

A



B



2.4. NEBNext End Prep

2.4.1. Use the Qubit readings from Step 2.3.13. to dilute 50 ng of the Targeted cDNA Amplicons sample with nuclease-free water to a final volume of 12.5 μl (4 ng/ μl). Add the following components to a PCR tube (NEBNext Ultra II End Prep Reaction Buffer and NEBNext Ultra II End Prep Enzyme Mix can be pre-mixed and this master mix is stable on ice for 4 hours):

COMPONENT	VOLUME
Targeted cDNA Amplicons (2.3.12)	12.5 μl
● (green) NEBNext Ultra II End Prep Reaction Buffer	1.75 μl
● (green) NEBNext Ultra II End Prep Enzyme Mix	0.75 μl
Total Volume	15 μl

2.4.2 Flick the tube or pipet up and down 10 times to mix the solution. 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 performance.

- 2.4.3 Place in a thermal cycler, with the heated lid set to $\geq 75^{\circ}\text{C}$, and run the following program:
 10 minutes @ 20°C
 10 minutes @ 65°C
 Hold at 4°C



If necessary, samples can be stored at -20°C for a few days; however, a slight loss in yield ($\sim 20\%$) may be observed. We recommend continuing with barcode ligation before stopping.

2.5. Barcode Ligation

- 2.5.1. Add the following components directly to a sterile nuclease-free PCR tube:

COMPONENT	VOLUME
○ (white) Nuclease-free water	6 μl
End-prepped DNA (2.4.3.)	1.5 μl
Native Barcode*	2.5 μl
● (red) Blunt/TA Ligase Master Mix**	10 μl
Total Volume	20 μl

* Native Barcodes are provided in Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104) and 13-24 (EXP-NBD114) or 1-96 (EXP-NBD196)

** Mix the Blunt/TA Ligase Master Mix by pipetting up and down several times prior to adding to the reaction

- 2.5.2. Flick the tube or pipet up and down 10 times to mix solution. Perform a quick spin to collect all liquid from the sides of the tube. **(Caution: The Blunt/TA Ligase Master Mix is very viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**
- 2.5.3. Incubate at room temperature for 20 minutes.
 Incubate at 65°C for 10 minutes.
 Place on ice for 1 min.
- 2.5.4. Pool all barcoded samples into one 1.5 ml DNA LoBind Tube.

2.6. Cleanup of Barcoded DNA

The following section is for cleanup of the ligation reaction.

Note: The volumes of SPRIselect or NEBNext Sample Purification Beads provided here are for use with the sample contained in the exact buffer at this step. AMPure XP Beads can be used as well. If using AMPure XP Beads, allow the beads to warm to room temperature for at least 30 minutes before use.

- 2.6.1. Vortex NEBNext Sample Purification Beads to resuspend.
- 2.6.2. Add 0.4X resuspended beads to pooled barcoded samples (2.5.4.) (for example, if you are pooling 24 barcoded DNA samples [which amounts to 480 μl total], add 192 μl of resuspended sample purification beads to the 480 μl of pooled sample). Flick the tube or pipet up and down 10 times to mix. Perform a quick spin for 1 second to collect all liquid from the sides of the tube.
- 2.6.3. Incubate samples on bench top for 10 minutes at room temperature.
- 2.6.4. Place the tube on a 1.5 ml magnetic stand (such as NEB #S1506) to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.
- 2.6.5. After 2 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets **(Caution: do not discard beads)**.
- 2.6.6. Wash the beads by adding 250 μl of Short Fragment buffer (SFB). Flick the tube or pipet up and down 10 times to mix to resuspend pellet. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 2.6.7. Place the tube on an appropriate magnetic stand for 2 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.

- 2.6.8. Repeat Step 2.6.6. and 2.6.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube, place back on the magnetic stand and remove traces of SFB with a p10 pipette tip
- 2.6.9. Add 200 μ l of 80% freshly prepared ethanol to the tube while on the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets.
- 2.6.10. A quick spin and place the sample tube on the magnetic stand, remove any residual ethanol.
- 2.6.11. Air dry the beads for 30 seconds 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. Elute the samples when the beads are still dark brown and glossy looking, but when all visible liquid has evaporated. When the beads turn lighter brown and start to crack, they are too dry.

- 2.6.12. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 33 μ l of Nuclease-free water.
- 2.6.13. Resuspend the pellet by flicking the tube or pipet up and down 10 times to mix. Incubate for at least 2 minutes at room temperature. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube before placing back on the magnetic stand.
- 2.6.14. Place the tube on the magnetic stand. After 2 minutes (or when the solution is clear), transfer 32 μ l to a new 1.5 ml Eppendorf DNA LoBind Tube or PCR tube.
- 2.6.15. Assess the concentration of the purified barcoded DNA sample. We recommend using a Qubit fluorometer for concentration assessment. Nanodrop is NOT recommended since it may overestimate the DNA concentration. Use 1 μ l of sample for the Qubit fluorometer.



Samples can be stored at -20°C if they are not used immediately.

2.7. Adapter Ligation

- 2.7.1. Use the Qubit readings from Step 2.6.15. to dilute up to 60 ng of the Native barcoded DNA pool with nuclease-free water to a final volume of 30 μ l (2 ng/ μ l). Add the following components into a 1.5 ml Eppendorf DNA LoBind Tube or nuclease-free PCR tube:

COMPONENT	VOLUME
Native barcoded and purified DNA (2.6.14.; up to 60 ng)	30 μ l
Adapter Mix II (AMII)**	5 μ l
• (red) NEBNext Quick Ligation Reaction Buffer*	10 μ l
• (red) NEBNext Quick T4 Ligase	5 μ l
Total Volume	50 μ l

* Mix the NEBNext Quick Ligation Reaction Buffer by pipetting up and down several times prior to adding to the reaction.

** Adapter Mix II is provided by Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104), 13-24 (EXP-NBD114) and 1-96 (EXP-NBD-196) kits.

- 2.7.2. Flick the tube to mix solution. Perform a quick spin for 1 second to collect all liquid from the sides of the tube. **(Caution: The NEBNext Quick Ligation Buffer is viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**
- 2.7.3. Incubate at 25°C or room temperature for 20 minutes.
- 2.7.4. Proceed to Cleanup of Adapter-ligated DNA in Section 2.8.

2.8. Cleanup of Adapter Ligated DNA

Note: The volumes of SPRIselect or NEBNext Sample Purification Beads provided here are for use with the sample contained in the exact buffer at this step. AMPure XP beads can be used as well. If using AMPure XP beads, allow the beads to warm to room temperature for at least 30 minutes before use. These volumes may not work properly for a cleanup at a different step in the workflow.

- 2.8.1. Vortex NEBNext Sample Purification Beads to resuspend.
- 2.8.2. Add 50 μ l (1X) resuspended beads to the ligation mix. Mix well by flicking the tube to mix followed by a quick spin for 1 second.

- 2.8.3. Incubate samples for 10 minutes at room temperature.
- 2.8.4. Place the tube on an appropriate magnetic stand to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.
- 2.8.5. After 2 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard the beads**).
- 2.8.6. Wash the beads by adding 250 μ l of Short Fragment Buffer (SFB). Flick the tube to resuspend pellet. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand. Place the tube on an appropriate magnetic stand.
- 2.8.7. Wait for 2 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.
- 2.8.8. Repeat steps 2.8.6. and 2.8.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube/plate, place back on the magnet and remove traces of SFB with a p10 pipette tip.
- 2.8.9. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 15 μ l of Elution Buffer (EB) provided in SQK-LSK114 kit from Oxford Nanopore.
- 2.8.10. Resuspend the pellet well in EB buffer by flicking the tube to mix. Incubate for 10 minutes at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 2.8.11. Place the tube/plate on the magnetic stand. After 2 minutes (or when the solution is clear), transfer 15 μ l to a new DNA LoBind tube.
- 2.8.12. Use Qubit to quantify 1 μ l DNA sample. Follow Oxford Nanopore Protocol SQK-LSK114 to prepare MinION[®] flow cell and DNA library sequencing mix using up to 20 ng adapter-ligated cDNA sample (2.8.11).

Note: After normalizing the DNA to 20 ng, if the volume is less than 12 μ l, then top off the sample volume to 12 μ l with EB.

Chapter 3

NEBNext VarSkip Short v2 Express Protocol without PCR Bead Cleanup (Two clean-up steps)

Symbols



This is a point where you can safely stop the protocol.



This caution sign signifies a step in the protocol that has two paths leading to the same end point but is dependent on a user variable, like the amount of input DNA.



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

Note 1: The amount of RNA required for detection depends on the abundance of the RNA of interest. In general, we recommend, using > 10 copies of the SARS-CoV-2 viral genome as input. The presence of carry-over products can interfere with sequencing accuracy, particularly for low copy targets. Therefore, it is important to carry out the appropriate no template control (NTC) reactions to demonstrate that positive reactions are meaningful. We recommend that the NTC reactions are prepared prior to non-NTC samples following the guidelines outlined below.

Guidelines for Sample Setup: We recommend setting up the cDNA synthesis and cDNA PCR setup reactions in a room (and ideally in a hood) separate from cDNA amplification and library construction area to minimize cross-contamination of future reactions.

Note 2: If sample Ct is between 12–15, then it is recommended per the ARTIC network *nCoV 2019 sequencing protocol v3 LoCost* to dilute the sample 100-fold in water, if between 15–18 then dilute 10-fold in water. This will reduce the likelihood of PCR-inhibition.

3.1. cDNA Synthesis

3.1.1. Gently mix and spin down the LunaScript RT SuperMix reagent. Prepare the cDNA synthesis reaction as described below:

COMPONENT	VOLUME
RNA Sample	8 μ l
● (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

3.1.2. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

3.1.3. For no template controls, mix the following components:

COMPONENT	VOLUME
○ (white) Nuclease-free Water	8 μ l
● (lilac) LunaScript RT SuperMix	2 μ l
Total Volume	10 μ l

3.1.4. Flick the tube or pipet up and down 10 times to mix followed by a quick spin.

3.1.5. Incubate reactions in a thermal cycler* with the following steps:

CYCLE STEP	TEMP	TIME
Primer Annealing	25°C	2 minutes
cDNA Synthesis	55°C	20 minutes
Heat Inactivation	95°C	1 minute
Hold	4°C	∞

*Set heated lid to 105°C



Samples can be stored at -20°C for up to a week.

3.2. Targeted cDNA Amplification

Note: 4.5 µl cDNA input is recommended. If using less than 4.5 µl of cDNA, add nuclease-free water to a final volume of 4.5 µl. NEBNext VarSkip Short v2 Primer Mixes cannot be combined with NEBNext ARTIC Primer Mixes in the same targeted amplification reaction.

Use of the NEBNext ARTIC Human Control Primer Pairs 1 and 2 are optional. If used, the appropriate NEBNext ARTIC Human Control Primer Pairs and NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix should be combined prior to use. More specifically, NEBNext ARTIC Human Control Primer Pairs 1 should be combined with NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1 and NEBNext ARTIC Human Control Primer Pairs 2 with NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2. Mixing directions are listed below.

3.2.1. Gently mix and spin down reagents. Prepare the split pool cDNA amplification reactions as described below:

For Pool Set A:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 1 with 42 µl of NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1 in a new tube, vortex and spin down reagents. If using a 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Mix 1 with 168 µl of NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1 in a new tube, vortex and spin down reagents. Use 1.75 µl of the combined mix for each Pool Set A reaction.

COMPONENT	VOLUME
cDNA (Step 3.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 1, add 1.75 µl of the combined VarSkip Short v2 SARS-CoV-2 Primer Mix 1 and NEBNext ARTIC Human Control Primer Pairs 1.

For Pool Set B:

If using the NEBNext ARTIC Human Control Primer Pairs and a 24 reaction kit, combine 0.7 µl of NEBNext ARTIC Human Control Primer Pairs 2 with 42 µl of NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2 in a new tube, vortex and spin down reagents. If using a 96 reaction kit, combine 2.8 µl of the NEBNext ARTIC Human Control Primer Pairs 2 with 168 µl of VarSkip Short SARS-CoV-2 Primer Mix 2 in a new tube, vortex and spin down reagents. Use 1.75 µl of the combined mix for each Pool Set B reaction.

COMPONENT	VOLUME
cDNA (Step 3.1.5)	4.5 µl
● (lilac) Q5 Hot Start High-Fidelity 2X Master Mix	6.25 µl
NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2*	1.75 µl
Total Volume	12.5 µl

* If using NEBNext ARTIC Human Control Primer Pairs 2, add 1.75 µl of the combined VarSkip Short v2 SARS-CoV-2 Primer Mix 2 and NEBNext ARTIC Human Control Primer Pairs 2.

Note: We recommend proceeding with the cDNA amplification and library construction in a different area or room

3.2.2. Incubate reactions in a thermal cycler* with the following steps:

CYCLE STEP	TEMP	TIME	CYCLES
Initial Denaturation	98°C	30 seconds	1
Denature	95°C	15 seconds	35
Annealing/Extension	63°C	5 minutes	
Hold	4°C	∞	1

3.2.3. Combine the Pool A and Pool B PCR reactions for each sample.



3.2.4. Samples can be stored at -20°C for up to a week.

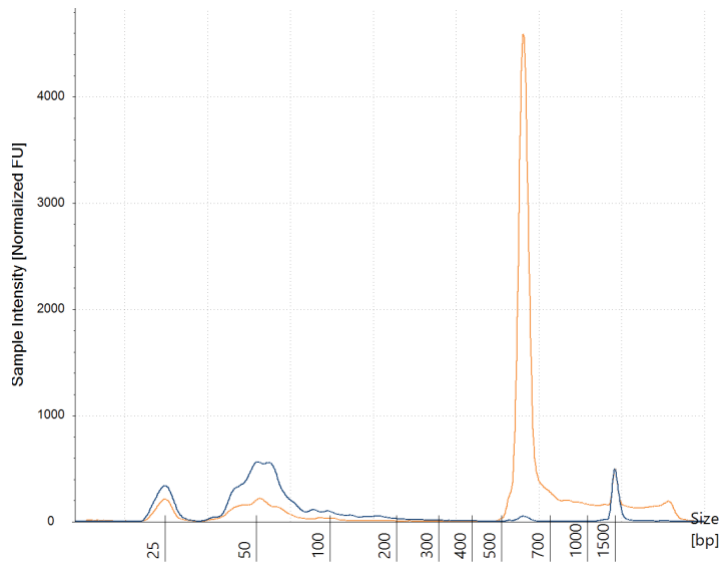
3.3. PCR Reaction Pooling

Please note, performing the cleanup and normalization step creates library pools where the reads for each library are more evenly distributed. These pools will likely achieve sufficient coverage in less run time

3.3.1. For each sample, combine pool A and pool B PCR Reactions.



Figure 3.3.1: VarSkip Short v2 SARS-CoV-2 cDNA amplicons generated from 1000 total viral copies. 1/10 diluted cDNA amplicons without bead cleanup run on a TapeStation in orange, NTC in blue.



3.4. NEBNext End Prep

3.4.1. Add the following components to a PCR tube (End Prep Reaction and Buffer can be pre-mixed and master mix is stable on ice for 4 hours):

COMPONENT	VOLUME
Targeted cDNA Amplicons (3.3.1)	1 µl
○ (white) Nuclease-free water	11.5 µl
● (green) NEBNext Ultra II End Prep Reaction Buffer	1.75 µl
● (green) NEBNext Ultra II End Prep Enzyme Mix	0.75 µl
Total Volume	15 µl

3.4.2. Flick the tube or pipet up and down 10 times to mix the solution. 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 performance.

3.4.3. Place in a thermal cycler, with the heated lid set to $\geq 75^{\circ}\text{C}$, and run the following program:

10 minutes @ 20°C

10 minutes @ 65°C

Hold at 4°C



If necessary, samples can be stored at -20°C for a few days; however, a slight loss in yield (~20%) may be observed. We recommend continuing with barcode ligation before stopping.

3.5. Barcode Ligation

3.5.1. Add the following components directly to a sterile nuclease-free PCR tube:

COMPONENT	VOLUME
○ (white) Nuclease-free water	6 μ l
End-prepped DNA (3.4.3.)	1.5 μ l
Native Barcode*	2.5 μ l
● (red) Blunt/TA Ligase Master Mix**	10 μ l
Total Volume	20 μ l

* Native Barcodes are provided in Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104) and 13-24 (EXP-NBD114) or 1-96 (EXP-NBD196)

** Mix the Blunt/TA Ligase Master Mix by pipetting up and down several times prior to adding to the reaction

3.5.2. Flick the tube or pipet up and down 10 times to mix solution. Perform a quick spin to collect all liquid from the sides of the tube. **(Caution: The Blunt/TA Ligase Master Mix is very viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**

3.5.3. Incubate at 25°C for 20 minutes.

Incubate at 65°C for 10 minutes.

Place on ice for 1 min.

3.5.4. Pool all barcoded samples into one 1.5 ml DNA LoBind Tube.

3.6. Cleanup of Barcoded DNA

Note: Use the pooled barcoded DNA samples (from Step 3.5.4.), up to 480 μ l for bead cleanup. Remaining pooled DNA can be stored at -20°C

3.6.1. Vortex NEBNext Sample Purification Beads to resuspend.

3.6.2. Add 0.4X resuspended beads to pooled barcoded samples (3.5.4.) (for example, if you are pooling 24 barcoded DNA samples [which amounts to 480 μ l total], add 192 μ l of resuspended sample purification beads to the 480 μ l of pooled sample). Mix well by flicking the tube or pipetting up and down 10 times. Perform a quick spin for 1 second to collect all liquid from the sides of the tube.

3.6.3. Incubate samples on bench top for 10 minutes at room temperature.

3.6.4. Place the tube on a 1.5 ml magnetic stand (such as NEB #S1506) to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.

3.6.5. After 3 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets **(Caution: do not discard beads).**

3.6.6. Wash the beads by adding 250 μ l of Short Fragment buffer (SFB). Flick the tube or pipet up and down 10 times to mix to resuspend pellet. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.

3.6.7. Place the tube on an appropriate magnetic stand for 3 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.

3.6.8. Repeat steps 3.6.6. and 3.6.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube, place back on the magnetic stand and remove traces of SFB with a p10 pipette tip

3.6.9. Add 500 μ l of 80% freshly prepared ethanol to the tube while on the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets. Repeat the wash once to make it a total of two washes.

3.6.10. A quick spin and place the sample tube on the magnetic stand, remove any residual ethanol.

3.6.11. Air dry the beads for 30 seconds 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. Elute the samples when the beads are still dark brown and glossy looking, but when all visible liquid has evaporated. When the beads turn lighter brown and start to crack, they are too dry.

3.6.12. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 33 μ l of Nuclease-free water.

- 3.6.13. Resuspend the pellet by flicking the tube or pipetting up and down 10 times to mix. Incubate for at least 2 minutes at room temperature. If necessary, quickly spin the sample for 1 second to collect the liquid from the sides of the tube before placing back on the magnetic stand.
- 3.6.14. Place the tube on the magnetic stand. After 2 minutes (or when the solution is clear), transfer 32 μl to a new 1.5 ml Eppendorf DNA LoBind Tube or PCR tube.
- 3.6.15. Assess the concentration of the purified barcoded DNA sample. We recommend using a Qubit fluorometer for concentration assessment. Nanodrop is NOT recommended since it may overestimate the DNA concentration. Use 1 μl of sample for the Qubit fluorometer.



Samples can be stored at -20°C if they are not used immediately.

3.7. Adapter Ligation

- 3.7.1. Use the Qubit readings from Step 3.6.15. to dilute up to 90 ng of the Native barcoded DNA pool with nuclease-free water to a final volume of 30 μl (3 ng/ μl). Add the following components into a 1.5 ml Eppendorf DNA LoBind Tube or nuclease-free PCR tube:

COMPONENT	VOLUME
Native barcoded and purified DNA (3.6.14.; up to 90 ng)	30 μl
• (red) NEBNext Quick Ligation Reaction Buffer*	10 μl
Adapter Mix II (AMII)**	5 μl
• (red) NEBNext Quick T4 Ligase	5 μl
Total Volume	50 μl

* Mix the NEBNext Quick Ligation Reaction Buffer by pipetting up and down several times prior to adding to the reaction.

** Adapter Mix II is provided by Oxford Nanopore Technologies Native Barcoding Expansion 1-12 (EXP-NBD104), 13-24 (EXP-NBD114) and 1-96 (EXP-NBD-196) kits.

- 3.7.2. Flick the tube to mix solution. Perform a quick spin for 1 second to collect all liquid from the sides of the tube. (**Caution: The NEBNext Quick Ligation Buffer is viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance.**)
- 3.7.3. Incubate at 25°C or room temperature for 20 minutes.
- 3.7.4. Proceed to Cleanup of Adapter-ligated DNA in Section 3.8.

3.8. Cleanup of Adapter Ligated DNA

Note: The volumes of SPRIselect or NEBNext Sample Purification Beads provided here are for use with the sample contained in the exact buffer at this step. AMPure XP beads can be used as well. If using AMPure XP beads, allow the beads to warm to room temperature for at least 30 minutes before use. These volumes may not work properly for a cleanup at a different step in the workflow.

- 3.8.1. Vortex NEBNext Sample Purification Beads to resuspend.
- 3.8.2. Add 50 μl (1X) resuspended beads to the ligation mix. Mix well by flicking the tube followed by a quick spin for 1 second.
- 3.8.3. Incubate samples for 10 minutes at room temperature.
- 3.8.4. Place the tube on an appropriate magnetic stand to separate the beads from the supernatant. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing on the magnetic stand.
- 3.8.5. After 3 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard the beads.**)
- 3.8.6. Wash the beads by adding 250 μl of Short Fragment Buffer (SFB). Flick the tube to resuspend pellet. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand. Place the tube on an appropriate magnetic stand.
- 3.8.7. Wait for 3 minutes (or until the solution is clear) to separate the beads from the supernatant. Remove the supernatant.

- 3.8.8. Repeat steps 3.8.6. and 3.8.7. once for a total of two washes. Be sure to remove all visible liquid after the second wash. If necessary, briefly spin the tube/plate, place back on the magnet and remove traces of SFB with a p10 pipette tip.
- 3.8.9. Remove the tube from the magnetic stand. Elute the DNA target from the beads by adding 15 μ l of Elution Buffer (EB) provided in SQK-LSK114 kit from Oxford Nanopore.
- 3.8.10. Resuspend the pellet well in EB buffer by flicking. Incubate for 10 minutes at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube or plate wells before placing back on the magnetic stand.
- 3.8.11. Place the tube/plate on the magnetic stand. After 3 minutes (or when the solution is clear), transfer 15 μ l to a new DNA LoBind tube.
- 3.8.12. Use Qubit to quantify 1 μ l DNA sample. Follow Oxford Nanopore Protocol SQK-LSK114 to prepare MinION[®] flow cell and DNA library sequencing mix using up to 30 ng adapter-ligated DNA sample (Step 3.8.11).

Note: After normalizing the DNA to 30 ng, if the volume is less than 12 μ l, then top off the sample volume to 12 μ l with EB.

Kit Components

NEB #E7660S Table of Components

NEB #	PRODUCT	VOLUME
E7651A	LunaScript RT SuperMix	0.048 ml
E7652A	Q5 Hot Start High-Fidelity 2X Master Mix	0.30 ml
E7661A	NEBNext Ultra II End Prep Enzyme Mix	0.018 ml
E7662A	NEBNext Ultra II End Prep Reaction Buffer	0.042 ml
E7663A	Blunt/TA Ligase Master Mix	0.24 ml
E7664A	NEBNext Quick T4 DNA Ligase	0.020 ml
E7665A	NEBNext Quick Ligation Reaction Buffer	0.040 ml
E7725A	NEBNext ARTIC SARS-CoV-2 Primer Mix 1	0.042 ml
E7726A	NEBNext ARTIC SARS-CoV-2 Primer Mix 2	0.042 ml
E7727A	NEBNext ARTIC Human Control Primer Pairs 1	0.007 ml
E7728A	NEBNext ARTIC Human Control Primer Pairs 2	0.007 ml
E7667A	Nuclease free-Water	1.50 ml
E7666S	NEBNext Sample Purification Beads	0.872 ml
E7872A	NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1	0.042 ml
E7872A	NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2	0.042 ml

NEB #E7660L Table of Components

NEB #	PRODUCT	VOLUME
E7651AA	LunaScript RT SuperMix	0.192 ml
E7652AA	Q5 Hot Start High-Fidelity 2X Master Mix	1.2 ml
E7661AA	NEBNext Ultra II End Prep Enzyme Mix	0.072 ml
E7662AA	NEBNext Ultra II End Prep Reaction Buffer	0.168 ml
E7663AA	Blunt/TA Ligase Master Mix	0.96 ml
E7664A	NEBNext Quick T4 DNA Ligase	0.020 ml
E7665A	NEBNext Quick Ligation Reaction Buffer	0.040 ml
E7725AA	NEBNext ARTIC SARS-CoV-2 Primer Mix 1	0.168 ml
E7726AA	NEBNext ARTIC SARS-CoV-2 Primer Mix 2	0.168 ml
E7727A	NEBNext ARTIC Human Control Primer Pairs 1	0.007 ml
E7728A	NEBNext ARTIC Human Control Primer Pairs 2	0.007 ml
E7667AA	Nuclease free-Water	4.7 ml
E7666L	NEBNext Sample Purification Beads	2.90 ml
E7872AA	NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1	0.168 ml
E7872AA	NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 2	0.168 ml

NEBNext ARTIC Human Primers

PRIMER MIX	GENE	POSITION	PRIMERS
NEBNext ARTIC Human Control Primer Pairs 1	EDF1	113 bp – 501 bp	GGCCAAATCCAAGCAGGCTA GTGTTTCATTTCCGCCCTAGGC
NEBNext ARTIC Human Control Primer Pairs 2	NEDD8	110 bp – 489 bp	AAAGTGAAGACGCTGACCGG GGGATCCTCACAGTCTCCCA

Detailed information for the ARTIC Human control primers can be found at: <https://doi.org/10.5281/zenodo.4495958>

NEBNext ARTIC SARS-CoV-2 Primer Mix 1 and 2

NEBNext ARTIC SARS-CoV-2 Primers for SARS-CoV-2 genome amplification are based on hCoV-2019/nCoV-2019 Version 3 (v3) sequences with balanced primer concentrations. Sequence information can be found at:

https://github.com/joshquick/artic-ncov2019/blob/master/primer_schemes/nCoV-2019/V3/nCoV-2019.tsv

NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1 and 2

NEBNext VarSkip Short v2 SARS-CoV-2 Mix 1 and 2 for SARS-CoV-2 genome amplification were designed to reduce the impact of variants on amplification efficiency. Sequence information can be found at:

<https://github.com/nebiolabs/VarSkip>

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	2/21
2.0	Update protocol	3/21
3.0	Update protocol and add second protocol	4/21
4.0	Update protocols	5/21
5.0	Add Chapter 3. Update protocols	9/21
6.0	Replace NEBNext VarSkip Short SARS-CoV-2 Primer Mix 1 and 2 for NEBNext VarSkip Short v2 SARS-CoV-2 Primer Mix 1 and 2.	2/22
7.0	Product change to the kit format.	3/22
8.0	Replace the discontinued ONT product SQK-LSK109 with the new version SQK-LSK114. Update format	5/26

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.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed. See www.neb.com/trademarks. The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email busdev@neb.com.

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. For additional information or to inquire about commercial use, please contact busdev@neb.com.

OXFORD NANOPORE TECHNOLOGIES® and MINION® are registered trademarks of Oxford Nanopore Technologies, LLC

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

AGILENT® and BIOANALYZER® are registered trademarks of Agilent Technologies, Inc.

AMPURE® is a registered trademark of Beckman Coulter, Inc.

B CORPORATION® is a registered trademark of B Lab IP, LLC, Inc.

THERMO FISHER SCIENTIFIC® is a registered trademark of Thermo Fisher Scientific Inc.

ALPAQUA® is a registered trademark of Alpaqua Engineering, LLC

EPPENDORF® is a registered trademark of Eppendorf AG

© Copyright 2026, New England Biolabs, Inc.; all rights reserved

