

Automated Preparation of Ion AmpliSeq™ SARS-CoV-2 Insight Research Panel Libraries

Tecan™ Fluent™ 1080 Automation Workstation

Publication Number MAN0025705 Revision B.0

■ Description	2
■ Laser safety scenarios	2
■ Required materials	5
■ Recommended materials	7
■ Before first use—Import script	7
■ Sample preparation guidelines	7
■ Prepare Instrument	9
■ Start the Tecan Fluent	11
■ Reverse transcribe RNA with the Ion Torrent™ NGS Reverse Transcription Kit	11
■ Prepare cDNA target amplification reactions	12
■ Partially digest amplicons	13
■ Ligate adapters	14
■ Purify the unamplified library	15
■ Equalize the library	15
■ Store libraries	16
■ Guidelines for templating and sequencing	17
■ Supplemental information	17
■ Documentation and support	19

IMPORTANT! This user bulletin is designed for experienced users of the Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay and Tecan™ Fluent™ 1080 Automation Workstation. For additional information, see the *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915) and *FluentControl Manual* (Pub. No. BG/N 30135092.04).

Note: For safety and biohazard guidelines, see the “Safety” appendix in the following product documentation: *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915) and *FluentControl Manual* (Pub. No. BG/N 30135092.04). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

This user bulletin describes how to prepare Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay – GS Manual libraries using Tecan™ Fluent™ 1080 Automation Workstation. The workflow for library preparation described in this bulletin is similar to the Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay workflow. Additional steps to set up the Tecan™ Fluent™ 1080 Automation Workstation and import and run the scripts are described.

For more information about the Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay, see the *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915).

For detailed instructions for using the Tecan™ Fluent™ 1080 Automation Workstation, see the *FluentControl Manual* (Pub. No. BG/N 30135092.04), available by contacting the Tecan Group at <https://lifesciences.tecan.com/>.

Laser safety scenarios



WARNING! LASER. Exposure to direct or reflected laser light can burn the retina and leave permanent blind spots. Never look into the laser beam. Remove jewelry and anything else that can reflect the beam into your eyes. Protect others from exposure to the beam.

For safety reasons, the lasers are disabled when they are outside of the CapSure™ LCM Cap area. However, there may be applications for which the cap is not required. In such instances, you can bypass the instrument laser safety settings.

To bypass the laser safety mechanism, depress the Laser Bypass button to activate the UV laser.

The following table details possible status scenarios, showing action combinations, and the resulting system responses. For example, row 4 indicates that when the laser bypass button has been pressed and a CapSure™ LCM Cap is in place, but the laser bypass key has *not* been inserted, the laser status is "Standby". You must insert the key into position for the laser to be ready to fire.

Table 1 Laser scenarios

State	Laser Bypass Button Pressed	Laser Bypass Key in Place	CapSure™ LCM Cap in Place	Laser State	To State	Pop-up	Cut Capture	Other	Tool Tip Messages
0 ^[1,2]				Standby			Stop		tt = 'Laser disabled because cap is out of beam path and override is off.'
1 ^[1, 2]				ON	0	x	Stop		tt = 'Laser disabled because cap is out of beam path and override is off. ' pop up = 'Lasers have been disabled because cap is out of beam path and override is off. Change cap placement or enable override to continue .'
2 ^[1,3]			x	Standby			Stop	Clean obj status	tt = 'Laser ready to fire. Cap placed in beam path.'
3 ^[4,5]			x	ON					tt = 'Laser is firing. Cap placed in beam path.'
4 ^[1,6]	x			Standby			Stop	Clean obj status	tt = 'Laser disabled because cap is out of beam path and override is off. '
5 ^[1,6]		x		ON	4	x	Stop	Clean obj status	tt = 'Laser disabled because cap is out of beam path and override is off. ' pop up = 'Lasers have been disabled because cap is out of beam path and override is off. Change cap placement or enable override to continue.'

Table 1 Laser scenarios *(continued)*

State	Laser Bypass Button Pressed	Laser Bypass Key in Place	CapSure™ LCM Cap in Place	Laser State	To State	Pop-up	Cut Capture	Other	Tool Tip Messages
6 ^[1,3]		x	x	Standby			Stop	Clean obj status	tt = 'Laser ready to fire. Cap placed in beam path.'
7 ^[4,5]		x	x	ON					tt = 'Laser is firing. Cap placed in beam path.'
8 ^[1,2]	x			Standby	0	x	Stop	Clean obj status	tt = 'Laser disabled because cap is out of beam path and override is off.' pop up = 'Lasers cannot be enabled because the hardware bypass key is absent.'
9 ^[1,2]	x			ON	0	x	Stop	Clean obj status	tt = 'Laser disabled because cap is out of beam path and override is off.' pop up = 'Lasers cannot be enabled because the hardware bypass key is absent.'
10									tt = 'Laser ready to fire. Cap placed in beam path.'
11 ^[4,5]	x		x	ON	3	x			tt = 'Laser is firing. Cap placed in beam path.' pop up = 'Laser interlock cannot be bypassed because the hardware key is not detected.'
12 ^[1,7]	x	x		Standby			Stop	Clean Obj status	tt = 'Laser ready to fire. Cap overridden and not placed in beam path.'
13 ^[4,8]	x	x		ON					tt = 'Laser is firing. Cap overridden and not detected in beam path.'

Table 1 Laser scenarios (continued)

State	Laser Bypass Button Pressed	Laser Bypass Key in Place	CapSure™ LCM Cap in Place	Laser State	To State	Pop-up	Cut Capture	Other	Tool Tip Messages
14 ^[1,9]	x	x	x	Standby			Stop	Clean obj status	tt = 'Laser ready to fire. Cap overridden but detected in beam path.'
15 ^[4,10]	x	x	x	ON					tt = 'Laser is firing. Cap overridden but detected in beam path.'

^[1] Set Laser: Standby

^[2] Color: U. Red

^[3] Color: U. Yellow

^[4] Set Laser: Power =p

^[5] Color: U. Green

^[6] Color: U. Orange

^[7] Color: PF. Yellow

^[8] Color: PF. Green

^[9] Color: P. Yellow

^[10] Color: P. Green

Required materials

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Item	Source
Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay – GS Manual	A51305
Ion Torrent™ NGS Reverse Transcription Kit	A45003
Instruments and Equipment	
Tecan™ Fluent™ 1080 Automation Workstation	http://tecan.com
One of the following thermal cyclers: <ul style="list-style-type: none"> GeneAmp™ PCR System 9700^[1] or GeneAmp™ PCR System 9700 96-Well^[1] 2720 Thermal Cycler^[1] Veriti™ 96-Well Thermal Cycler ProFlex™ 96-well PCR System 	See web product pages
MicroAmp™ Splash-Free 96-Well Base	4312063
MicroAmp™ Optical Film Compression Pad	4312639
Alpaqua™ 96S Super Magnet	A001322 (Alpaqua™)
96-well plate centrifuge	MLS

(continued)

Item	Source
Reagents and consumables	
150 µL MCA Disposable Tips	30180837 (Tecan)
MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode	4483354, 4483352
Thermo Scientific™ Nunc™ 96-Well Polypropylene DeepWell™ Storage Plates	12-565-395
MicroAmp™ Clear Adhesive Film	4306311
Agencourt™ AMPure™ XP Reagent	A63880, A63881, or A63882 (Beckman Coulter™)
Nuclease-free water	AM9932
70% v/v Ethanol solution	BP82011
Corning™ 96-well Clear V-Bottom 2 mL Polypropylene Deep Well Plate, Sterile	3960 (Corning™)
Agilent™ 300 mL Reservoir	201244100 (Agilent™)
One or more of the following kits for nucleic acid isolation and quantification	
RecoverAll™ Total Nucleic Acid Isolation Kit for FFPE	AM1975
MagMAX™ FFPE DNA/RNA Ultra Kit	A31881
PureLink™ Genomic DNA Mini Kit	K1820-00
(Recommended for DNA quantification) TaqMan™ RNase P Detection Reagents Kit	4316831
(Recommended for RNA quantification) Qubit™ RNA HS Assay Kit	Q32852 or Q32855
Barcodes	
Ion Xpress™ Barcode Adapters Kit	Various
IonCode™ Barcode Adapters 1–384 Kit	A29751
Ion Torrent™ Dual Barcode Kit 1–96	A39360
One or more of the following kits for quantification	
Ion Library TaqMan™ Quantitation Kit	4468802
Ion Library Equalizer™ Kit	4482298
If you are not using the Ion Library Equalizer™ Kit for library normalization, select one of the following kits:	
Qubit™ Fluorometer ^[2] and Qubit™ dsDNA HS Assay Kit	Q33238, Q32851 or Q32854
Agilent™ 2100 Bioanalyzer™ and Agilent™ High Sensitivity DNA Kit	G2939BA, 5067-4626 (Agilent™)

^[1] Supported but no longer available for purchase.

^[2] Qubit™ 2.0 Fluorometer or later

Recommended materials

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Item	Source
Recommended for SARS-CoV-2 Quantification	
TaqPath™ COVID-19 Combo Kit	A47814
Recommended for RNA dilution	
THE RNA Storage Solution	AM7000

Before first use—Import script

Scripts are instructions for the Tecan™ Fluent™ 1080 Automation Workstation. You must download the appropriate script for preparing Ion AmpliSeq™ SARS-CoV-2 Insight Research Panel libraries.

1. Open the Tecan™ FluentControl™ Software.
2. Select **Database ▶ Import....**
3. Select **AmpliSeq SARS-CoV-2 Insight Research Assay 780.zeia**.
4. Press **Add all ▶ Import**, remove any detected conflicts, then, if prompted, select **Yes** to import without the missing referenced files.

Sample preparation guidelines

Guidelines for sample quality, viral copy number, and variant calling

- The amount of viral RNA among samples should be approximately equivalent so that the target amplification conditions you select are optimal for all samples.
- Ensure that RNA samples are quantified using TaqPath™ COVID-19 Combo Kit (Cat. No. [A47814](#)).

- A sample containing as little as 50 copies of viral RNA after isolation (25 copies per target amplification reaction) can be used to prepare an Ion AmpliSeq™ SARS-CoV-2 Insight Research Panel library. For optimal results, we recommend a viral copy number in the 100 to 200,000 range, or an amount of *total* RNA between 1–10 ng.

Table 2 Sample quality and viral copy number

Viral copy number	Recommendations and guidelines
200 to 200,000 copies	Recommended range for optimal results.
50 to 200 copies	Only for high-quality samples without degradation. We recommend sequencing and variant detection with a minimum allele frequency of 20%. For more information about the minimum allele frequency, see the Torrent Suite™ Software Help.

- To reliably sequence low quality samples, the samples must have a viral copy number ≥ 200 copies per reaction. For partially degraded samples, which likely includes low titer samples, the effective copy number that can be amplified by the Ion AmpliSeq™ SARS-CoV-2 Insight Research Panel is lower than the viral copy number detected by qPCR because the qPCR products are shorter than the 250 bp fragments generated by the panel.
- Even for samples with viral titers >200 copies per reaction, you may observe reverse transcription-derived false positives if you decrease the minimum allele frequency cutoff below 0.2 (20%). Reverse transcription-related errors occur randomly across the genome. To minimize calling false-positives, be certain to amplify a sufficient number of RNA molecules and set the minimum allele frequency to at least 20%.
- See the *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915) for recommended RNA isolation and quantification kits.

Copy number determination by qPCR

- If your qPCR data give a different relationship between C_t and copy number, this is likely a result of differences in the baseline or threshold selected. Determine the copy number of a sample according to the known copy number in control reactions.
- We recommend basing copy number on the N Protein C_t value.
- If the N Protein C_t value is not accurate, use the S Protein or ORF1ab C_t values to determine copy number.
- The copy number is only an estimate.

Table 3 Approximate copy number to C_t conversion—TaqPath™ COVID-19 RT-PCR kits

Tier	Viral copy number	TaqPath™ C_t		
		N Protein	S Protein	ORF1ab
Low	50–1,500	25–29	24–29	24–29
Medium	1,500–50,000	20–25	19–24	19–24
High	50,000–1,500,000	15–20	14–19	15–19

Guidelines for RNA isolation, quantification, and input

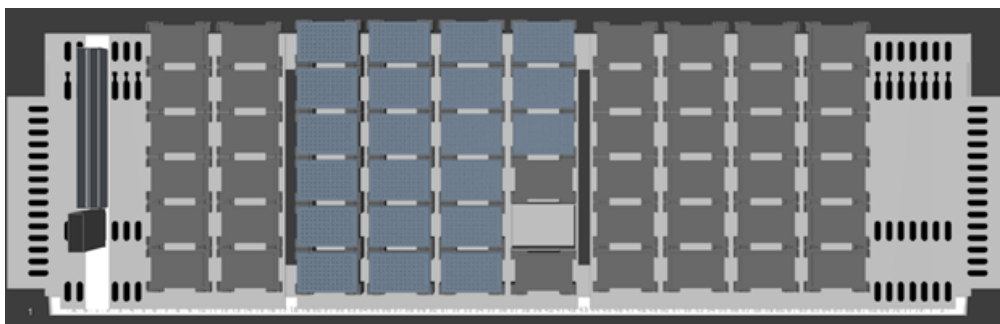
- For recommended kits for isolating RNA, see “Required materials” on page 5.
- Each reverse transcription reaction requires 1–100 ng of DNase-treated RNA (≥ 0.14 ng/ μ L), prepared from normal or formaldehyde- or paraformaldehyde-fixed paraffin-embedded (FFPE) tissue.
- For quantifying RNA, we recommend the Qubit™ RNA HS Assay Kit (Cat. No. [Q32852](#) or [Q32855](#)).
- In general, the library yield from high-quality RNA is greater than from degraded samples. Library yield is not indicative of sequencing performance.
- Having more RNA starting material generally results in higher quality libraries. However, if RNA is not degraded, high-quality libraries can be generated from as little as 1 ng starting material.

Prepare Instrument

Guidelines for Tecan Fluent

The procedure described in this user bulletin has many steps based on a specific configuration of a Tecan™ Fluent™ 1080 Automation Workstation. If you are using any other configuration you must ensure performance through simulated runs and wet runs using mock solutions before testing real samples.

Sets of sample plates are grouped according to standard Tecan nomenclature such that Labware ending in the same [xxx] are part of the same sample set. For example, RNA[001] is processed in RT[001] and RNA[002] is processed in RT[002]. The worktable has been setup to have sample groups placed within the same site position on neighboring grid segments.



Segment description	Grid location
Waste Thru Trough 8x100ml RL 00	1
(Optional) Segment Deck 3 Grids	3
6 Landscape 7mm Nest	6
6 Landscape 7mm Nest	12
Empty	18
6 Landscape 61mm Nest	19

(continued)

Segment description	Grid location
6 Landscape 61mm Nest	25
6 Landscape 61mm Nest	31
4 Landscape 61mm Nest Thru Deck Waste (with additional 61mm Nest at Site 1)	37
Empty	43
6 Landscape 7mm Nest	44
6 Landscape 7mm Nest	50
6 Landscape 7mm Nest	56
6 Landscape 7mm Nest	62

Prepare reagent master plates

Aliquot Ion AmpliSeq™ reagents into MicroAmp™ EnduraPlate™ Optical 96-Well Fast Clear Reaction Plates.

For details, see “Recommended fill volumes for reagent master plates (high throughput applications)” on page 18.

Recommended reagent substitutions for calibration

Use the following reagent substitutions during initial setup of the Tecan™ Fluent™ 1080 Automation Workstation for calibration. You can also use the recommended reagent substitutions for troubleshooting of any performance problems and inconsistencies.

Reagent	Recommended substitution
DNA	Water
Primers	
Ion Torrent™ NGS 10X RT Enzyme Mix	50% glycerol solution in water
Ion Torrent™ NGS 5X Reaction Buffer	
FuPa Reagent	
DNA Ligase	
5X Ion AmpliSeq™ HiFi Mix	40% glycerol solution in water
Switch Solution	20% Polyethylene Glycol 8000 (PEG-8000) solution in water
Agencourt™ AMPure™ XP Reagent	


Tip handling

The worktable is setup with 36 MCA tip boxes and is sufficient to complete all scripts without intervention. The MCA tips do not need to be replaced between runs. The script prompts if a tip refresh is required for the next run.

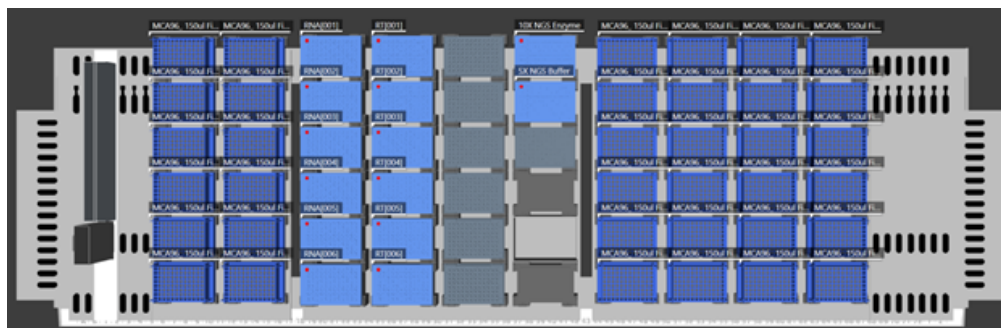
Start the Tecan Fluent

1. Power on the Tecan™ Fluent™ 1080 Automation Workstation.
2. Open Tecan™ FluentControl™ Software.
3. If prompted, enter your username and password.
4. Initialize the instrument by selecting **Run ▶ Initialize Instrument**.

Reverse transcribe RNA with the Ion Torrent™ NGS Reverse Transcription Kit

1. Follow the alternate reverse transcription protocol in the *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915).
2. In the Tecan™ FluentControl™ Software, touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Reverse Transcribe** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
3. Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.


The estimated run time for 6 plates is 11 minutes.



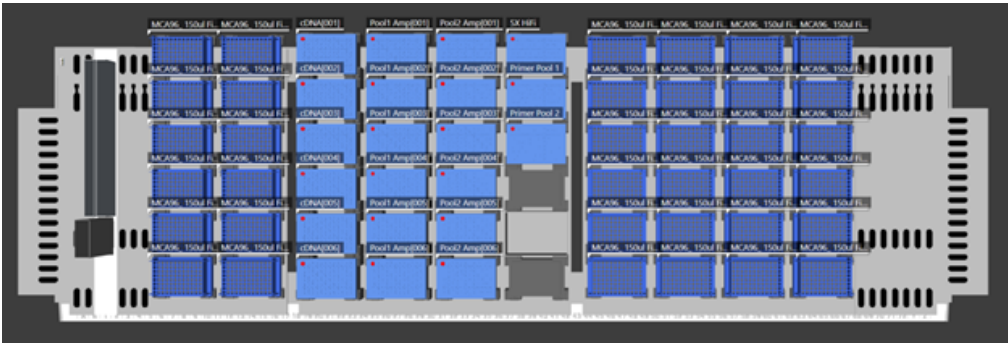
4. Run the following program to reverse transcribe RNA.

Temperature	Time
25°C	10 minutes
50°C	10 minutes
85°C	5 minutes
10°C	Hold

Prepare cDNA target amplification reactions

1. Touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Amplify cDNA** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
2. Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.

The estimated run time for 6 plates is 20 minutes.



3. Run the following program to amplify the target regions.

Stage	Step	Temperature	Time
Hold	Activate the enzyme	98°C	2 min
Cycle; set number according to Table 4	Denature	98°C	15 sec
	Anneal and extend	60°C	4 min
Hold	—	10°C	Hold

Table 4 Recommended cycle number

Tier	Viral copy number	Number of amplification cycles
Low	50–1,500	26
Medium	1,500–50,000	20
High ^[1]	50,000–1,500,000	15

^[1] If titers are above 1,500,000 copies, samples can be diluted.


Cycle number recommendations in the preceding table are based on qPCR quantification of viral copy number. Without qPCR quantification, use the following guidelines to determine optimal cycle number empirically.

- Low viral load suspected: 26 cycles.
- High viral load suspected: 20 cycles.
- Isolates or enriched viral particles: ~15 cycles for 2 ng input.

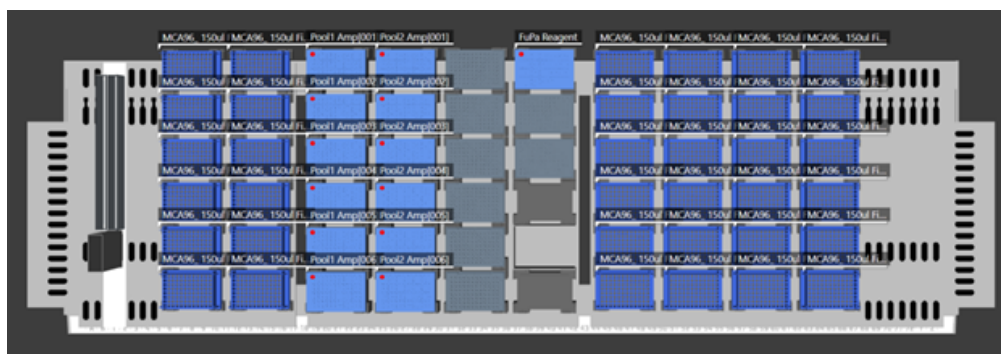
If you are working with samples with **unknown viral load**, and cannot quantify using qPCR, use 20 target amplification cycles as a starting point for manual library preparation.

STOPPING POINT Target amplification reactions can be stored at 10°C overnight on the thermal cycler. For longer periods, store at –20°C.

Partially digest amplicons

1. Touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Partially Digest** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
2. Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.


The estimated run time for 6 plates is 10 minutes.



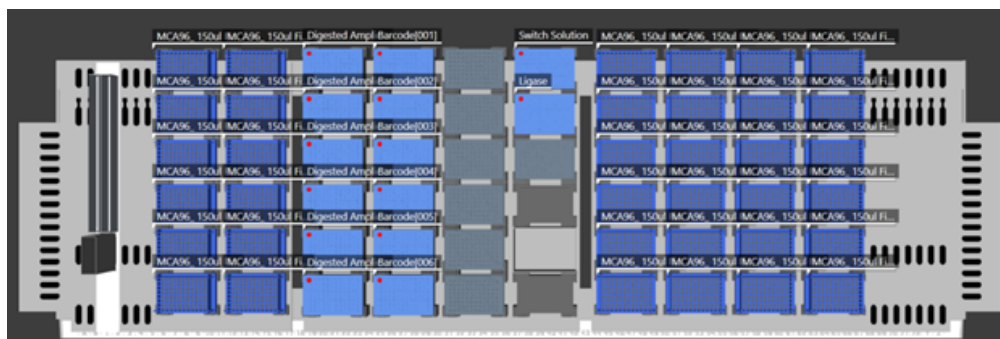
- Run the following program to amplify the target regions.

Temperature	Time
50°C	10 min
55°C	10 min
60°C	20 min
10°C	Hold (for up to 1 hour)

Ligate adapters

- Touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Ligate Adapters** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
- Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.


The estimated run time for 6 plates is 12 minutes.



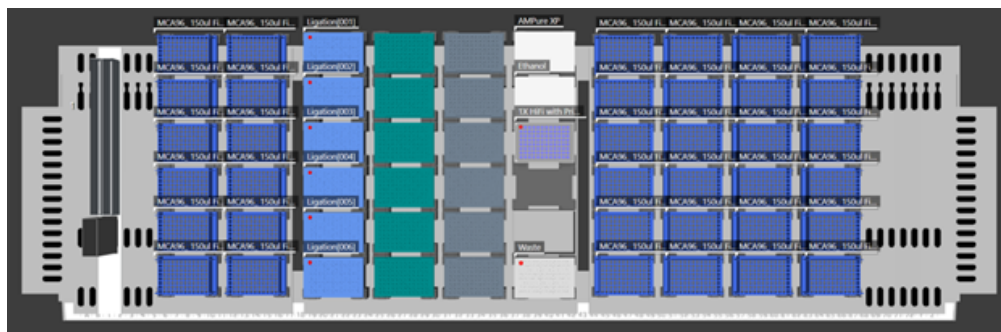
- Run the following program to amplify ligate adapters.

Temperature	Time
22°C	30 minutes
68°C	5 minutes
72°C	5 minutes
10°C	Hold (for up to 24 hours)

Purify the unamplified library

1. Touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Purify & Amplify** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
2. Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.


The estimated run time for 6 plates is 48 minutes.

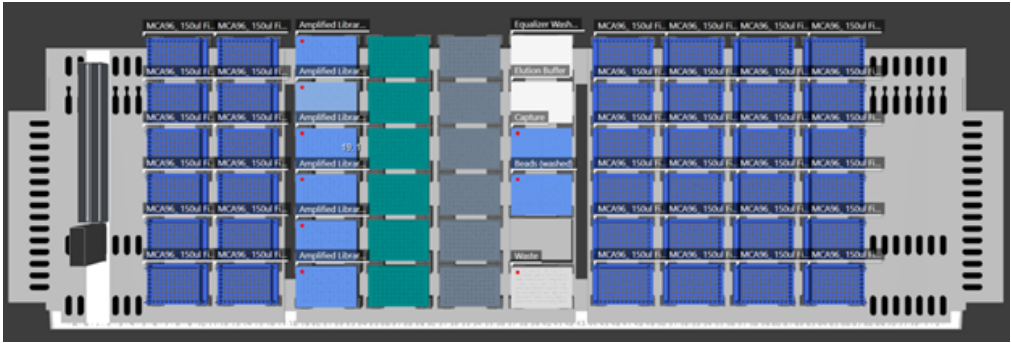


3. Run the following program to amplify the target regions.

Stage	Temperature	Time
Hold	98°C	2 minutes
9 cycles	98°C	15 seconds
	64°C	1 minute
Hold	10°C	Hold (up to 1 hour)

Equalize the library

1. Ensure that the Equalizer™ Beads have been washed. For instructions, see *Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide* (Pub. No. MAN0024915).
 2. Touch  (**Method Starter**), then select **AmpliSeq SARS-CoV-2 Equalize Library** to start the run. Alternatively, the method can be started by selecting the method within Tecan™ FluentControl™ Software.
 3. Using the touch screen, select the number of plates to run. Follow the prompts for placement of the plates and minimum required fill volumes for reagents. The setup and volumes are matched to the number of plates being processed.
- The estimated run time for 6 plates is 55 minutes.



4. Perform thermal cycling using the conditions in the following table.

Stage	Temperature	Time
Hold	32°C	5 minutes
Hold	10°C	Hold (up to 1 hour)

The supernatant contains the Equalized library at ~100 pM, which can be stored with beads for up to 1 month at 4–8°C.

5. Dilute library to the appropriate concentration.

Chip	Concentration
Ion 530™ Chip	30 pM
Ion 540™ Chip	50 pM

Proceed to templating and sequencing.

Store libraries

Libraries may be stored at 4–8°C for up to 1 month. For longer term storage, store at –20°C.

Guidelines for templating and sequencing

Proceed to template preparation and sequencing using the following kits.

Chip	Maximum libraries/chip	Kit	User Guide
Ion 530™ Chip ^[1,2]	<ul style="list-style-type: none"> 16^[3] 32^[4] 	Ion 510™ & Ion 520™ & Ion 530™ Kit – Chef (Cat. No. A34461)	<i>Ion 510™ & Ion 520™ & Ion 530™ Kit – Chef User Guide</i> (Pub. No. MAN0016854)
Ion 540™ Chip ^[1,2]	<ul style="list-style-type: none"> 64^[3] 128^[4] 	Ion 540™ Kit – Chef (Cat. No. A30011)	<i>Ion 540™ Kit – Chef User Guide</i> (Pub. No. MAN0010851)

^[1] Template system: Ion Chef™ System

^[2] Sequencer: Ion S5™ XL Sequencer, Ion GeneStudio™ S5 Plus Sequencer, or Ion GeneStudio™ S5 Prime Sequencer

^[3] 1,000,000 reads

^[4] 500,000 reads

Supplemental information

Minimum fill volumes required for each run configuration

The script prompts these fill values when starting a run. The following table shows the required volume per well (µL). Where possible, fill master plates with 12 reactions per well to minimize dead volume losses (see “Recommended fill volumes for reagent master plates (high throughput applications)” on page 18).

Component	Number of plates					
	1	2	3	4	5	6
Reverse transcription reagents						
Ion Torrent™ NGS 5X Reaction Buffer	8	11	14	17	20	23
Ion Torrent™ NGS 10X RT Enzyme Mix	6.5	8	9.5	11	12.5	14
cDNA target amplification reagents						
5X Ion AmpliSeq™ HiFi Mix	9	13	17	21	25	29
Ion AmpliSeq™ 5X primer pool	7	9	11	13	15	17
Amplicon digestion reagents						
FuPa Reagent	7	9	11	13	15	17
Ligation reagents						
DNA Ligase	7	9	11	13	15	17
Switch Solution	9	13	17	21	25	29

(continued)

Component	Number of plates					
	1	2	3	4	5	6
IonCode™, Ion Xpress™, or Ion Torrent™ Dual Barcode Adapters	6	6	6	6	6	6
Library purification reagents						
AMPure™ beads	18.5	22	25.5	29	32.5	36
70% Ethanol	45	75	105	135	165	195
Equalization reagents						
1X Ion AmpliSeq™ HiFi + primers	75	130	185	240	295	350
Equalizer™ Capture	15	25	35	45	55	65
Equalizer™ Beads, washed	11	17	23	29	35	41
Equalizer™ Wash Buffer	45	75	105	135	165	195
Equalizer™ Elution Buffer	25	35	45	55	65	75

Recommended fill volumes for reagent master plates (high throughput applications)

To minimize dead volume loss, we recommend filling reagent master plates to support up to 12 sample plates (1,152 total samples). If accessing the plates multiple times, ensure the total number of uses is tracked to avoid dry wells. The estimated uses per plate is 12.

Reagent	Dead volume (µL)	Recommended fill volume per well (µL)
5X Reaction Buffer	5	41
10X RT Enzyme Mix	5	23
5X Ion AmpliSeq™ HiFi Mix	5	53
Ion AmpliSeq™ SARS-CoV-2 Insight Research Panel Pool ^[1]	5	29
FuPa Reagent	5	29
DNA Ligase	5	29
Switch Solution	10	58
1X Library Amplification Mix + Equalizer primers ^[2]	20	680

(continued)

Reagent	Dead volume (µL)	Recommended fill volume per well (µL)
Equalizer™ Capture	5	125
Washed Equalizer™ Beads	5	71

[1] You must prepare a plate for each primer.

[2] Use the Thermo Scientific™ Nunc™ 96-Well Polypropylene DeepWell™ Storage Plates.

Documentation and support

Related documentation

Document	Publication number
<i>Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay User Guide</i>	MAN0024915
<i>Ion AmpliSeq™ Library Kit Plus User Guide</i>	MAN0017003
<i>FluentControl Manual</i>	BG/N 30135092.04 (https://lifesciences.tecan.com/)

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, California 92008 USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: MAN0025705 B.0 (English)

Revision	Date	Description
B.0	29 January 2024	<ul style="list-style-type: none">Removed references to discontinued TaqMan™ 2019-nCoV Assay Kit v1 (A47532).Removed references to discontinued TaqPath™ COVID-19 CE-IVD RT-PCR Kit (A51738/A48067).
A.0	18 February 2022	New user bulletin for preparing Ion AmpliSeq™ SARS-CoV-2 Insight Research Assay libraries on the Tecan™ Fluent™ 1080 Automation Workstation.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

TRADEMARKS: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tecan, Fluent, and FluentControl are trademarks of Tecan Group, Ltd. Bioanalyzer and Agilent are trademarks of Agilent Technologies, Inc. TaqMan is a trademark of Roche Molecular Systems, Inc., used under permission and license. Agencourt and AMPure are trademarks of Beckman Coulter, Inc. Alpaqua is a trademark of Alpaqua Engineering, LLC.

©2022-2024 Thermo Fisher Scientific Inc. All rights reserved.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

29 January 2024

ThermoFisher
SCIENTIFIC