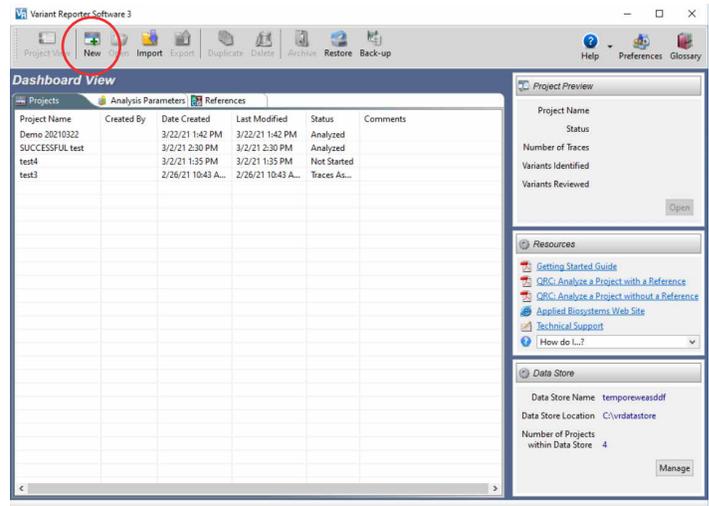


# Suggestions for SARS-CoV-2 variant analysis of Sanger sequencing traces using Variant Reporter Software

Detecting and confirming mutations in newly arising SARS-CoV-2 strains is an important strategy for controlling the spread of new viral lineages. Recently we published four protocols that use Sanger sequencing to analyze SARS-CoV-2 isolates (protocols available at [thermofisher.com/sangercoronavirus](https://thermofisher.com/sangercoronavirus)). Although the data contained in the sequence files can be compared against known sequences using a BLAST™ search, in some cases a simpler and more definitive workflow is needed. In this protocol, we present options for SARS-CoV-2 variant analysis using Sanger sequencing files and Applied Biosystems™ Variant Reporter™ Software. Although this workflow is not comprehensive, it is presented to help users begin to better understand their results. To take advantage of the complete features of the software, please see the user guide.

Setup of a basic project in Variant Reporter Software generally involves the following steps: importing sequencing traces (.ab1 files), defining a reference sequence, importing known variants of the reference sequence, and importing primer sequences defining the PCR amplicons.

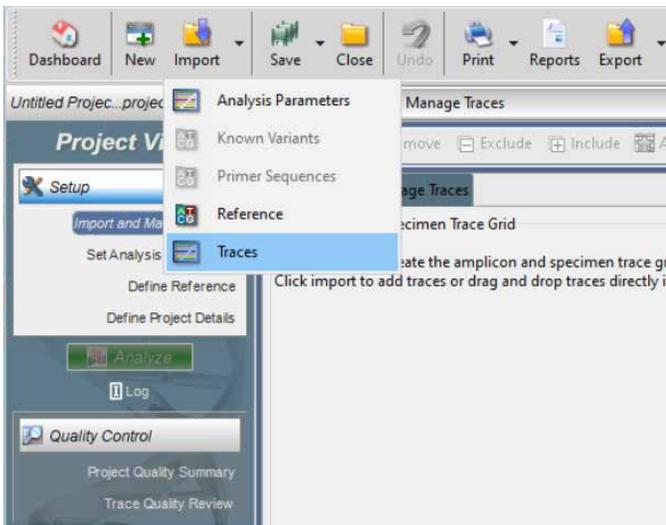
## 1. Create new project



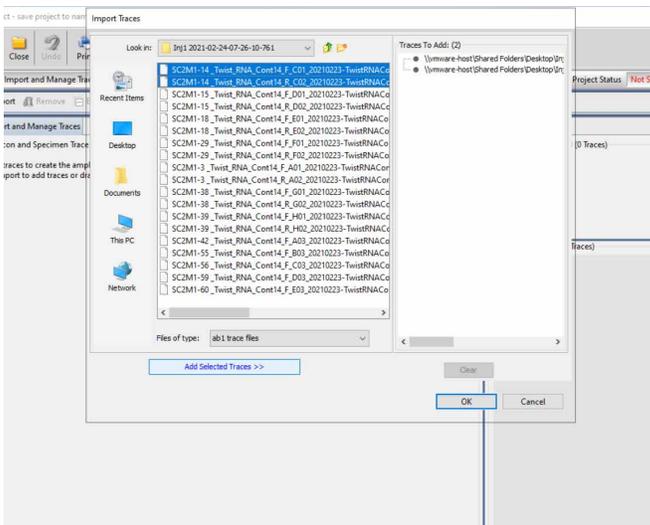
The Import and Manage Traces page is used to import traces into your project, creating the Amplicon and Specimen Trace Grid. From the application dashboard, click the New icon to create a new project.

## 2. Import sequencing traces

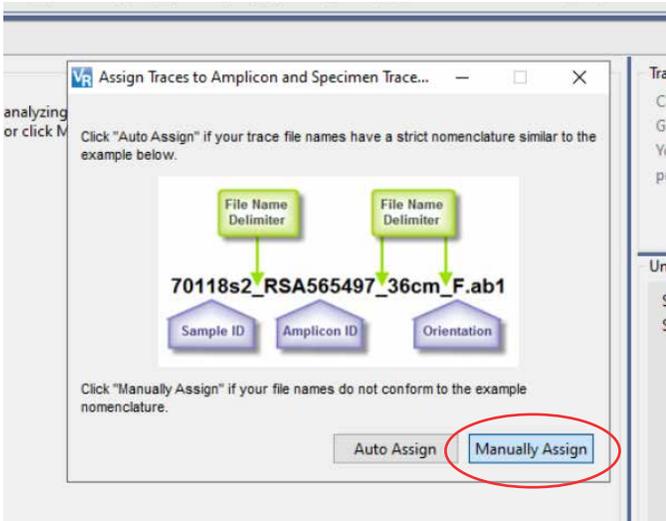
**A**



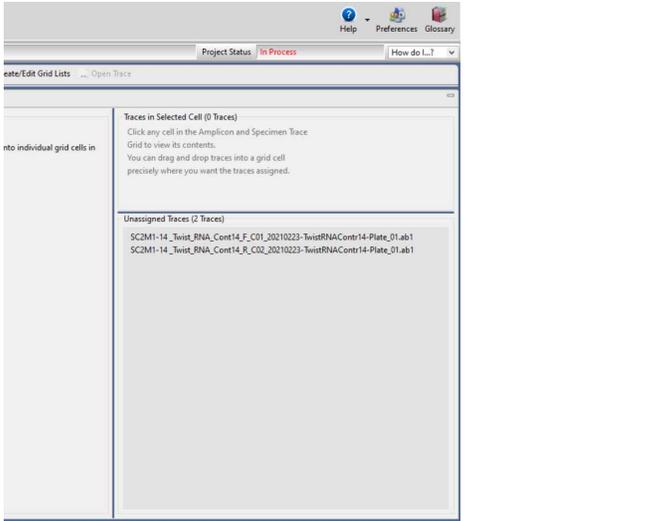
**B**



**C**



**D**



The figure consists of four panels (A, B, C, D) illustrating the process of importing sequencing traces into a software application.

**(A)** The software interface shows the 'Import' menu with 'Traces' selected. The 'Traces' option is highlighted in blue.

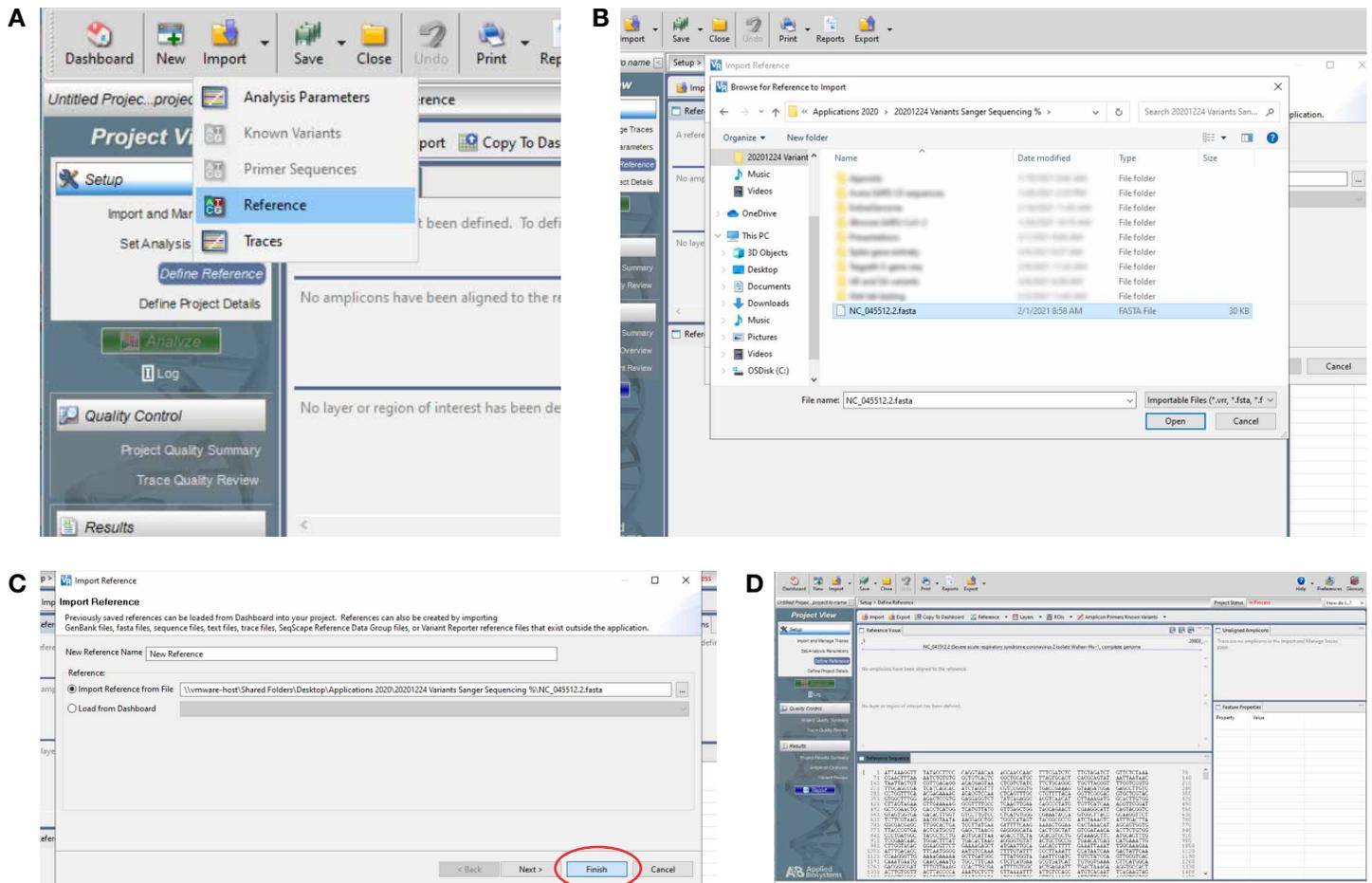
**(B)** The 'Import Traces' dialog box is shown, displaying a list of files to be imported. The files are listed in a table with columns for file name and file type. The file type is set to 'ab1 trace files'. The 'Add Selected Traces >>' button is highlighted.

**(C)** The 'Assign Traces to Amplicon and Specimen Trace...' dialog box is shown. It provides instructions on how to assign traces. An example file name '70118s2\_RSA565497\_36cm\_F.ab1' is shown, with arrows pointing to 'File Name Delimiter' and 'File Name Delimiter' labels. Below the example, three boxes represent 'Sample ID', 'Amplicon ID', and 'Orientation'. The 'Manually Assign' button is circled in red.

**(D)** The software interface shows the 'Unassigned Traces' section. The traces are listed in a table with columns for file name and file type. The file type is set to 'ab1 trace files'. The 'Unassigned Traces' section is highlighted.

**(A)** To import the sequencing traces, choose Traces from the Import menu. **(B)** Next, navigate to the folder containing the traces and choose all that are to be imported. Select all files to be analyzed, click Add Selected Traces, then click OK. **(C)** If the file name is in the appropriate format (see example above), the traces can be assigned sample names, amplicons, and sequencing direction automatically. However, for the purposes of this demonstration, the files will be assigned manually by clicking the Manually Assign button. **(D)** Once the button is clicked, traces that have been successfully imported but not yet assigned to amplicons will be listed in the bottom-right quadrant of the subsequent window.

### 3. Import reference sequence

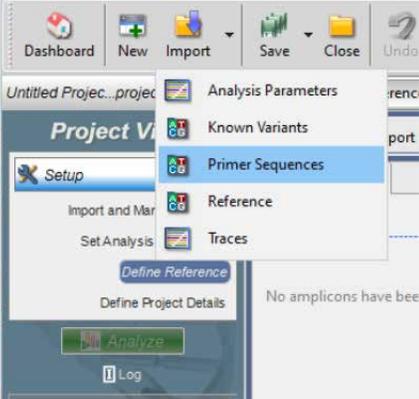


(A) To import the reference sequence, choose Reference from the Import menu. (B) Navigate to the reference text file. In this case, the SARS-CoV-2 reference strain NC\_045512.2 is used. Variant Reporter Software also supports reference sequences in the .gb file format to include sequence annotations. The FASTA file can be downloaded from [thermofisher.com/sangercoronavirus](https://thermofisher.com/sangercoronavirus). Choose the desired FASTA file and click Open. (C) A window appears confirming the location of the file and giving an opportunity to rename it. When completed, click Finish. (D) The imported sequence will appear in its own pane at the bottom of the window.

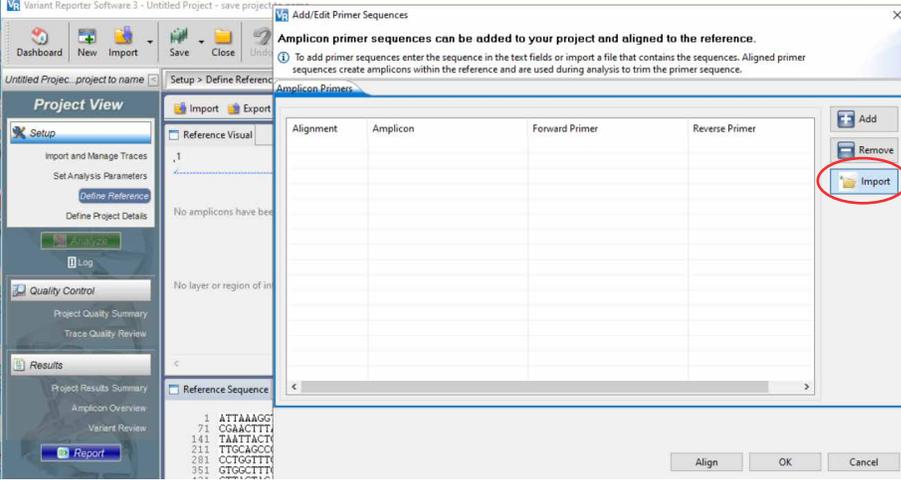


## 5. Import primer sequences that define amplicons used

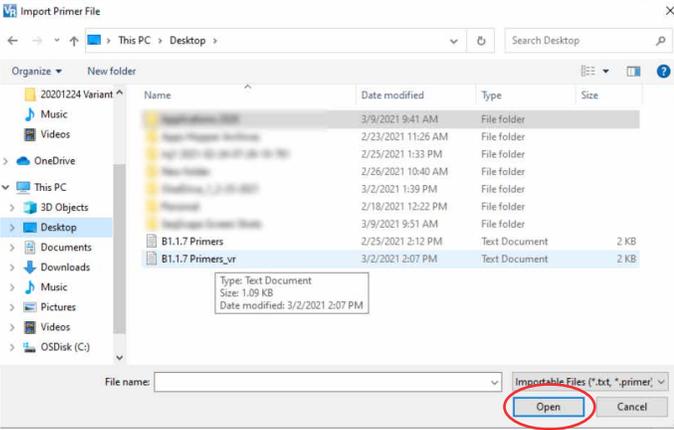
**A**



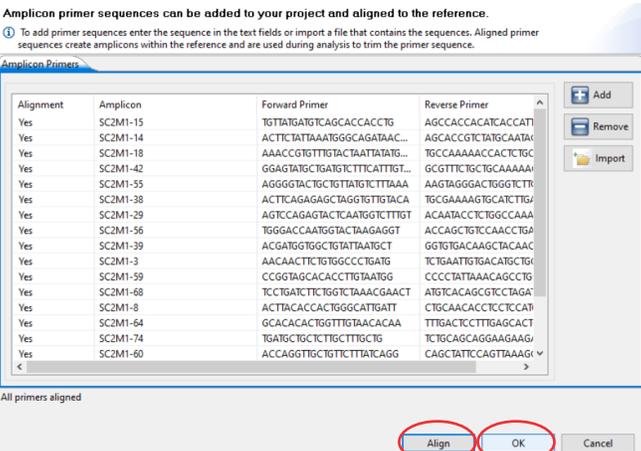
**B**



**C**



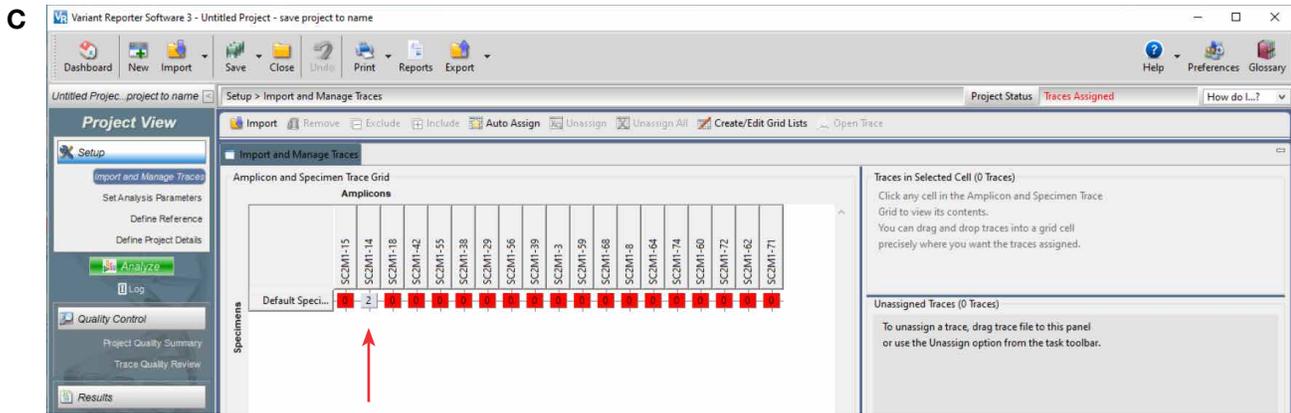
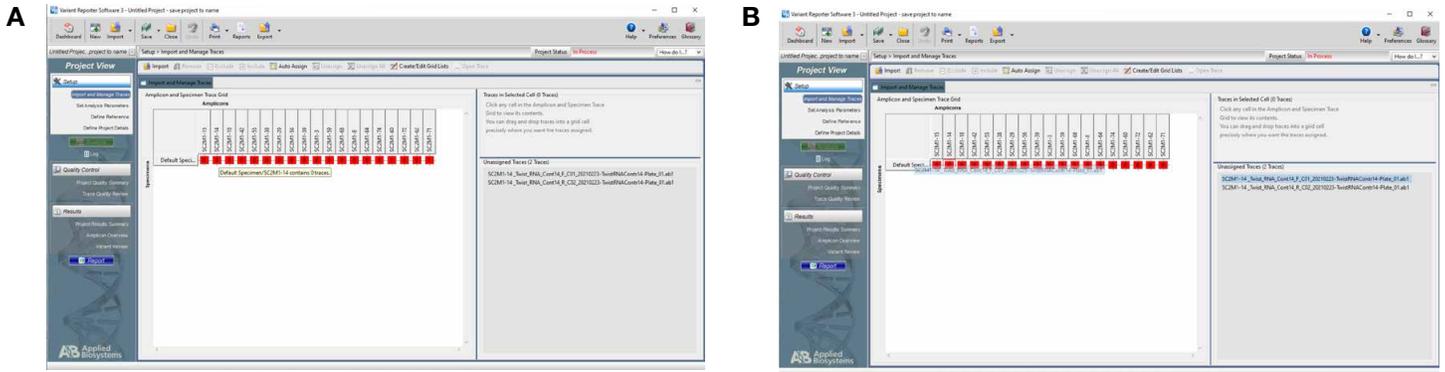
**D**



Alignment	Amplicon	Forward Primer	Reverse Primer
Yes	SC2M1-15	TGTTATGATGTCAGCCACCCTG	AGCCACCACATCACCATI
Yes	SC2M1-14	ACTTCTATTAATAATGGCAGATAAC...	AGCACCCTGTATGCAATAA
Yes	SC2M1-18	AAACCGTGTGTGACTAATATATG...	TGCCAAAAACCACTCTGC
Yes	SC2M1-42	GGAGTAGCTGATGTCCTTTCATTG...	GGCTTCTGCTGCAAAAAA
Yes	SC2M1-55	AGGGGTACTGCTGTATGTCCTTAA...	AAGTAGGGACTGGGCTCT
Yes	SC2M1-38	ACTTCAGAGAGCTGGGTTGTTTACA	TGCGAAAAGTGCATCTG
Yes	SC2M1-29	AGTCAGAGATGCTCAATGCTCTTGT	ACATATCTCTGCGCAAA
Yes	SC2M1-56	TGGACCAATGACTAAGAGGTT	ACCACCTGCTCAACCTG
Yes	SC2M1-39	ACCATGTTGGCTGATATGCT	GGTGTGCAAGTACAACT
Yes	SC2M1-3	AACAACCTCTGTGGCCCTGATG	TCGTAAATGTCATGCTG
Yes	SC2M1-59	CCGGTAGCACACCTGTAATGG	CCCTATTAACAGCCCTG
Yes	SC2M1-68	TCTGTATCTTGTCTAAACGAAC	ATGTACACAGGCTCTAG
Yes	SC2M1-8	ACTTACACCTGSGGATGAT	CTCACACCTCTCTCAT
Yes	SC2M1-64	GCACACACTGTTTAAACAGAA	TTTGACTCTTISAGACT
Yes	SC2M1-74	TGATGTCCTGCTGCTGCTG	TCGTGAGCAGGAAGAG
Yes	SC2M1-60	AACCGGTGCTGCTCTTATCAGG	CAGCTATCCAGTTAAAG

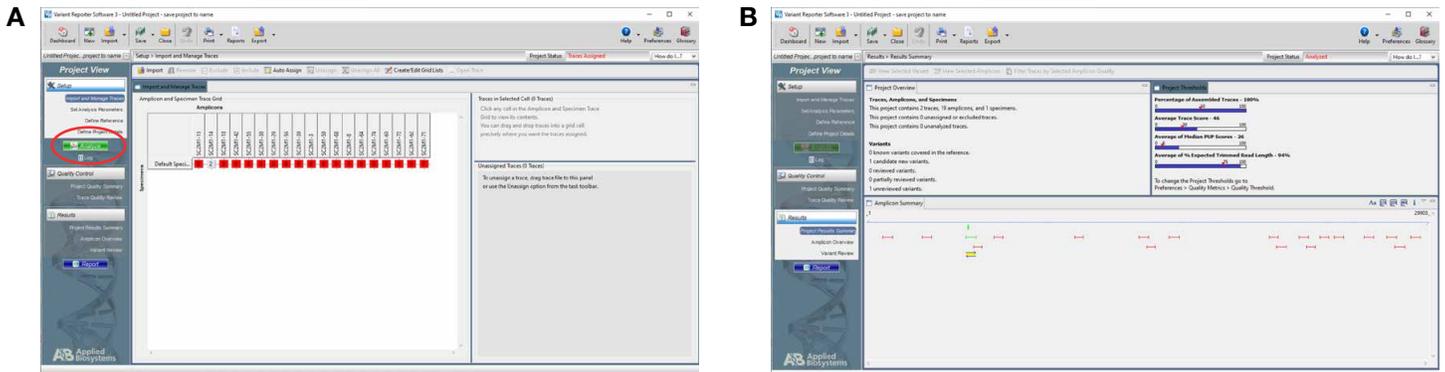
Next, the sequences of the primers defining the amplicons need to be imported. A text file containing primer sequences can be downloaded from [thermofisher.com/sangercoronavirus](https://thermofisher.com/sangercoronavirus). Download the file to a local computer. **(A)** From the Import menu, choose Primer Sequences. **(B)** A window appears; choose Import and navigate to the location of the downloaded text file. **(C)** Choose the file and click Open. **(D)** Once the primers have been imported, align them to the reference sequence by clicking the Align button. If the primers are found in the reference sequence, the alignment column will show Yes. When all primers have been imported successfully, click OK.

## 6. Assign traces to sequencing primers (manual)



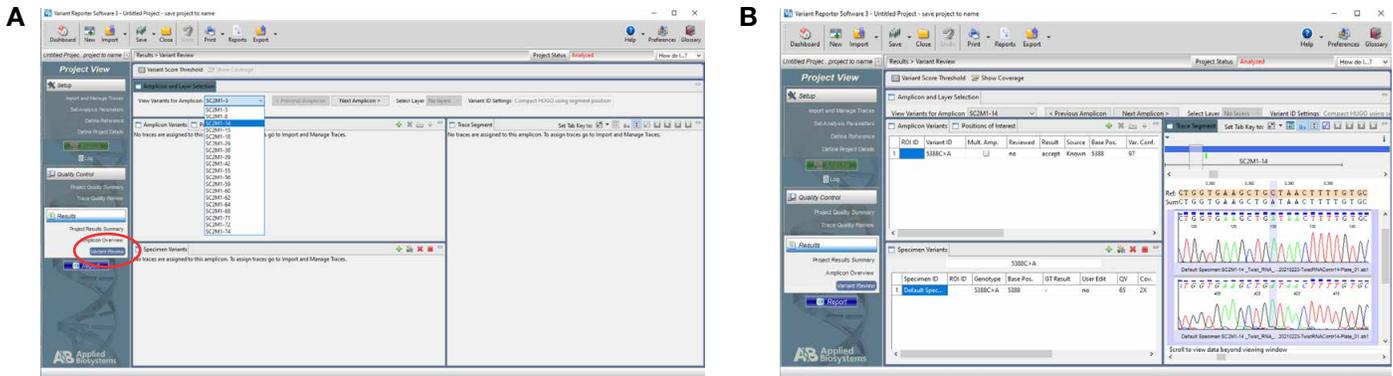
(A) When primers are imported, a grid showing all the primer pairs will appear. Next, sequencing traces will be assigned to the correct primer pairs in the grid. To assign the traces to the primers, drag and drop the file from the list on the right to the correct position in the grid on the left. (B) Here, we are moving the file SC2M1-14\_Twist\_RNA\_Cont14\_F\_C01\_20210223-TwistRNACont14-Plate\_01.ab1 to the square under the SC2M1-14 box in the Amplicon and Specimen Trace Grid. (C) The color of the box will change from red (empty) to yellow when one trace is added, then gray when the reverse trace is added. Forward and reverse traces from the same amplicon are added to the same square in the grid. Continue until all traces are assigned.

## 7. Analyze traces



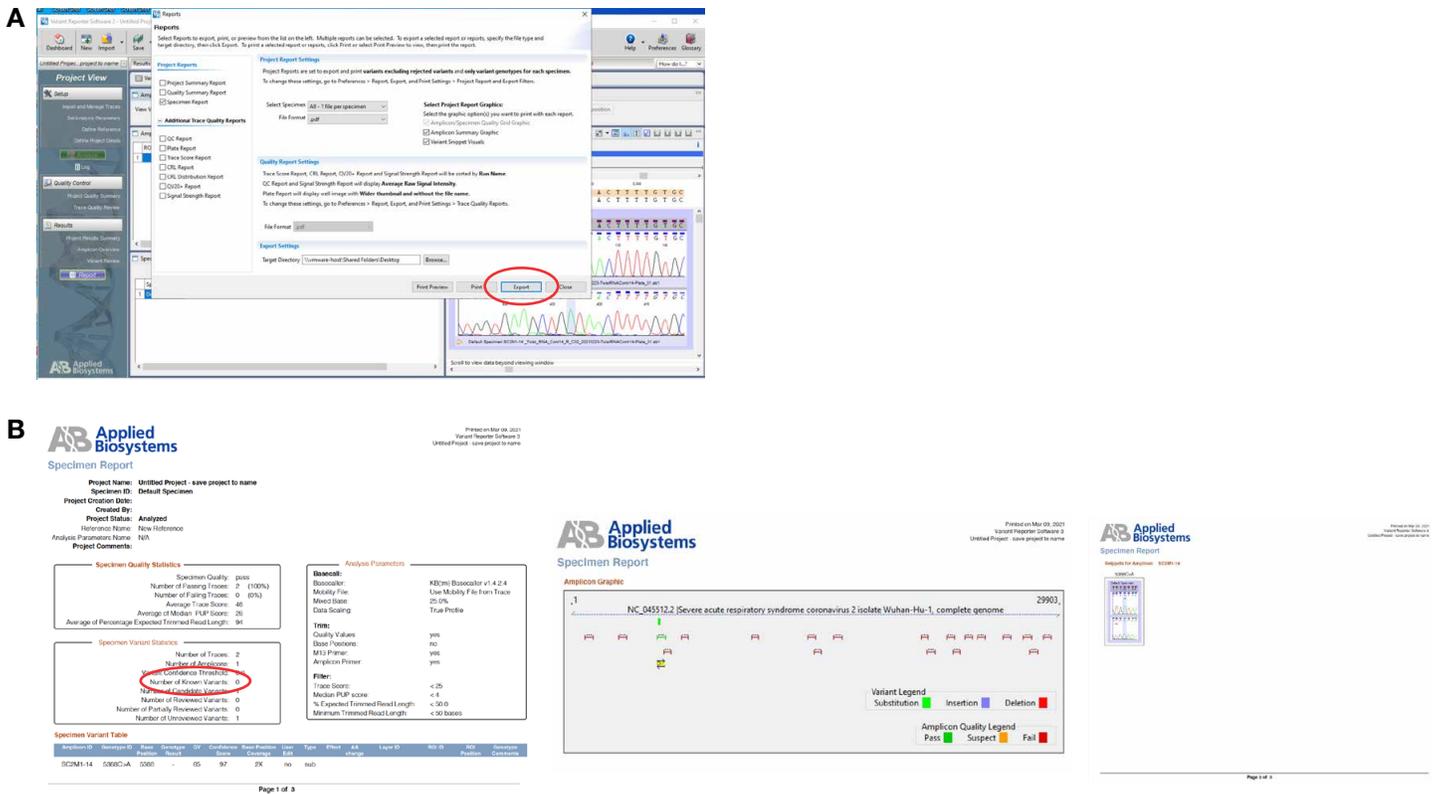
(A) To analyze the traces, click the green Analyze button on the left. (B) The position of the amplicon is shown in green, and the traces that were successfully analyzed are shown below the amplicon and highlighted in yellow. In addition, a summary of the analyzed traces is shown on the right side of the window.

## 8. Choose amplicon to view results



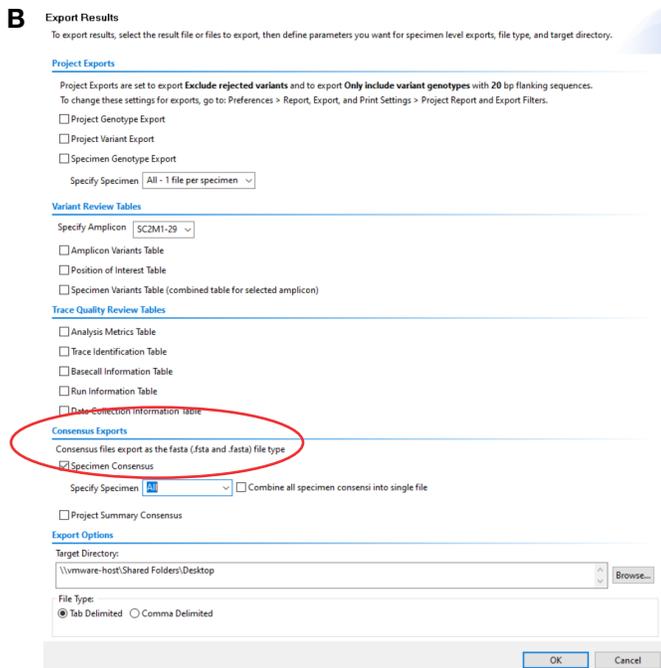
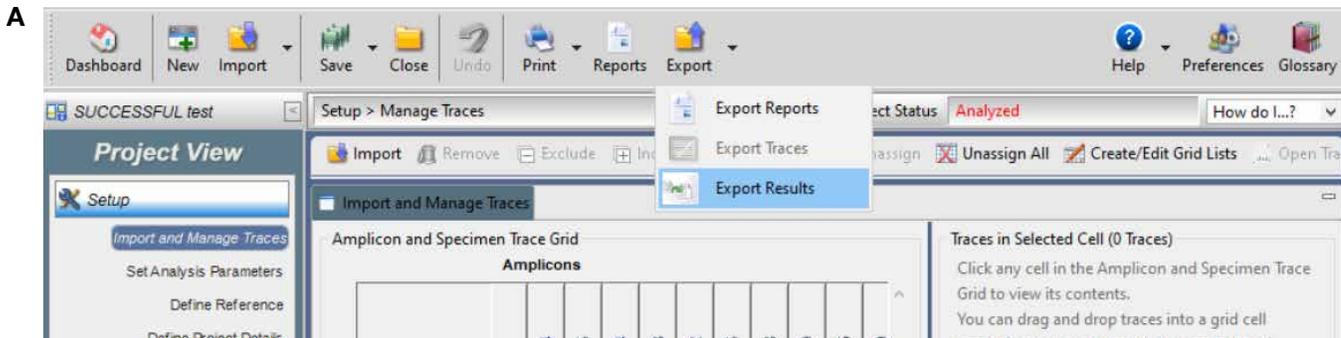
(A) Variants can be visualized by selecting Variant Review from the Project View panel on the left side of the screen (red circle). Then, select the amplicon from the drop-down menu as shown. (B) Variants that map to the amplicon will be shown in the next window. Here, the variants within the amplicon and in the sample are listed, as well as the electropherograms showing the sequence change (far right). Navigation between the variants can be done by clicking Next Amplicon at the top of the window.

## 9. Generate report



(A) A complete list of the results of the project can be generated by clicking the Reports menu on the main screen. A window with choices for the report appears. The report can be exported by choosing Specimen Report from the choices on the left side of the window, followed by Export. (B) Examples of the three pages from the resulting PDF are shown. If any known variants were found in the sample, they will appear on the first page of the report (circled).

## 10. Mapping sequences to known strain lineages



Results can be exported into FASTA files for querying public databases. **(A)** To generate one or more FASTA files, choose Export Results from the Export icon in the toolbar. **(B)** In the window that appears, choose Specimen Consensus near the bottom of the form (red circle). Choose which specimens to export, specify the directory where the exported files will appear, then click OK. A text file containing FASTA files for each of the amplicons in each specimen will be generated. Sequences in this file can be used to query strain lineage databases. One of these, the GISAID database ([epicov.org/epi3/frontend#637476](http://epicov.org/epi3/frontend#637476)), provides BLAST search results against all known SARS-CoV-2 strains. Up to 10 FASTA sequences can be queried at once. Registration is required, but use is free. The Coronavirus Typing Tool ([genomedetective.com/app/typingtool/cov](http://genomedetective.com/app/typingtool/cov)) also accepts FASTA files. Single FASTA files can be queried at no charge, but batch uploading requires premium access.

Find out more at [thermofisher.com/sangercoronavirus](http://thermofisher.com/sangercoronavirus)