

CarrierScan™ Reporter v1.1

USER GUIDE

Publication Number 703467

Revision 3



Affymetrix, Inc.
3450 Central Expressway
Santa Clara, CA 95051

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

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

Revision history: Pub No. 703467

| Revision | Date | Description |
|----------|--------------|---|
| 3 | June 2019 | Version 1.1 release with updated Licensing information. |
| 2 | May 2019 | Version 1.1 release |
| 1 | January 2018 | Initial release |

Important Software Licensing Information

Your installation and/or use of this CarrierScan Reporter software is subject to the terms and conditions contained in the End User License Agreement (EULA) which is incorporated within the CarrierScan Reporter software, and you will be bound by the EULA terms and conditions if you install and/or use the software.

Legal entity

Affymetrix, Inc. | Santa Clara, CA 95051 USA | Toll Free in USA 1 800 955 6288

TRADEMARKS

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

©2019 Thermo Fisher Scientific Inc. All rights reserved.

| | | |
|---|--|-----------|
| ■ | CHAPTER 1 Introduction | 5 |
| | Overview | 5 |
| | Features | 6 |
| | New features in v1.1 | 6 |
| | Minimum requirements | 6 |
| | Installation | 7 |
| | Start up | 7 |
| | Locating the library folder | 8 |
| | Locating the output folder | 8 |
| | (Optional) Creating copy number visualization files | 8 |
| | Setting up custom proxy settings | 8 |
| | Downloading library files | 8 |
| ■ | CHAPTER 2 Setup | 9 |
| | Importing data | 9 |
| | Selecting an analysis mode | 12 |
| | Manual mode | 12 |
| | Individual analysis (one sample) | 12 |
| | Adding multiple samples to an individual analysis simultaneously | 13 |
| | Paired analysis (two samples) | 14 |
| | Removing an analysis name from the setup | 14 |
| | Running an analysis | 14 |
| | Batch mode | 16 |
| | Required batch mode file guidelines | 16 |
| | Searching table content | 18 |
| | Removing all samples from the table | 19 |
| | Updating the library folder | 19 |
| ■ | CHAPTER 3 Summary tables | 20 |
| | Batch summary table | 20 |
| | Selecting previously generated results | 20 |
| | Displaying only pathogenic variants | 22 |
| | Summary table | 23 |
| | Searching table content | 24 |
| | Exporting the summary table | 24 |
| ■ | CHAPTER 4 Results table | 25 |
| | Using the results table | 25 |
| | Selecting a results folder | 25 |
| | Searching table content | 27 |
| | Default row sorting | 27 |

Contents

| | | |
|---|---|-----------|
| ■ | CHAPTER 5 Exporting | 28 |
| | Exports | 28 |
| | Full export | 28 |
| | Positive export | 30 |
| | Summary export | 30 |
| | Review export | 31 |
| ■ | APPENDIX A Chance of occurrence calculation guidelines | 32 |



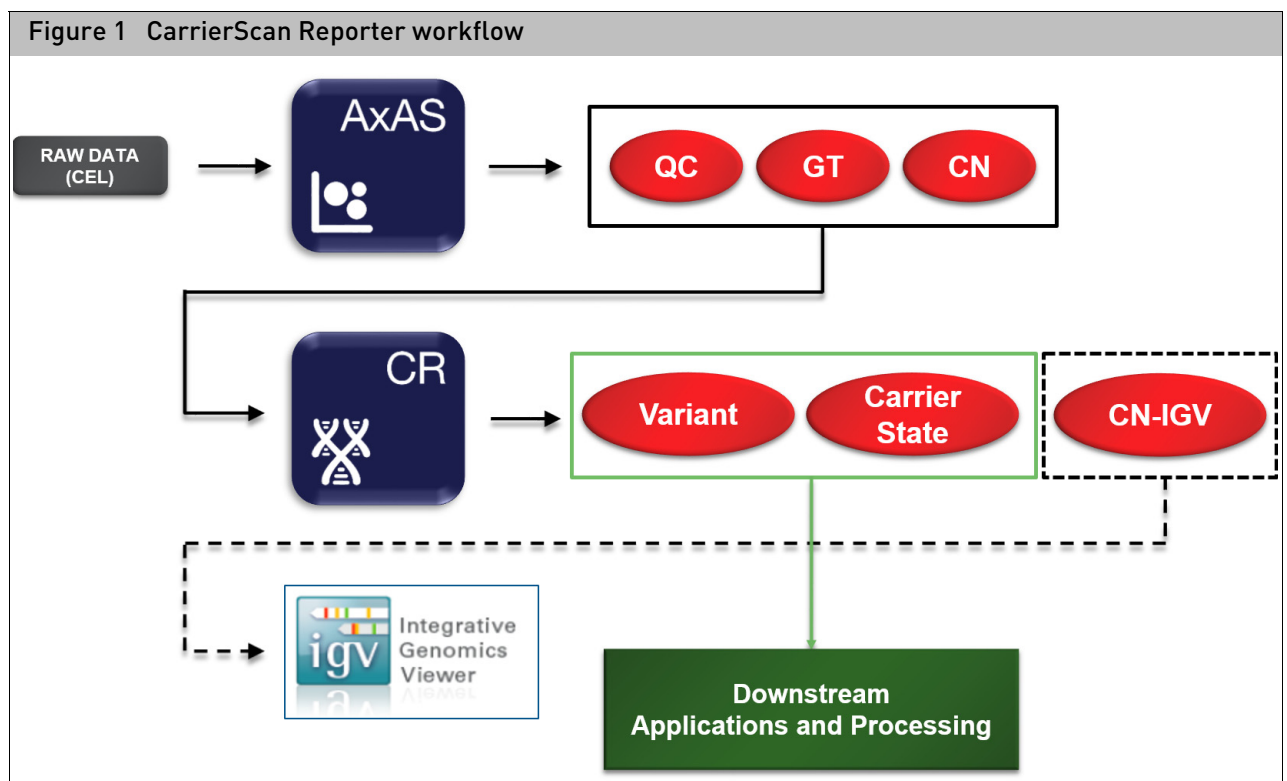
Introduction

Overview

CarrierScan Reporter is designed to work with Axiom Analysis Suite (AxAS) and its best practices analysis of CarrierScan based arrays. AxAS calculates quality control metrics, as well as assigns genotyping and copy number assessments. It reads in the AxAS results, then determines sample heredity status using phenotypic descriptions found in the public domain (ClinVar, OMIM, etc.).

CarrierScan Reporter also assesses individual and paired samples, then creates customizable reports of the annotated results.

The flowchart (Figure 1) describes the CarrierScan data generation pipeline from analyzing CEL files in Axiom Analysis Suite to CarrierScan Reporter variant and carrier translation. The text outputs can be imported and processed in existing analysis pipelines. In addition, all carrier determinations can be traced back to the original genotype and copy number calls and graphically assessed in AxAS and Integrative Genomics Viewer (IGV), respectively.



Features

- Controlled and automated sample carrier state determinations.
- Individual residual chance of carrier occurrence.
- Paired residual chance of phenotype occurrence.
- Compiled information from public databases for associated phenotypes.

New features in v1.1

- CarrierScan Reporter v1.1 uses a new annotation panel containing annotation content for the full array. Output will still be limited to the desired variants based on AxAS results. Custom annotation panels limiting to the desired content are no longer needed. **Note:** CarrierScan Reporter v1.1 requires new library files. See ["Downloading library files" on page 8](#).
- A sample QC metric "Rare Carrier Count" has replaced the QC Het Rate metric. This sample QC identifies samples with >20 rare variants called as non-normal and adjusts the calls to No Call as samples with a high number of rare carrier/affected calls is indicative of an issue with the sample.
- The requirement to have all annotation columns present in the annotation panel file has been removed. Any annotation column present in the annotation panel will be displayed, but is no longer required.
- Multi-allele variants are now exported on individual lines similar to bi-alleles for easier viewing.
- The link between HBA1 and HBA2 has been removed and will be exported individually.
- The output folder will automatically be named similar to the AxAS Batch name for easier matching of AxAS data and CarrierScan Reporter data for Carrier Reviewer Tool.

Minimum requirements

| Operating System | Processor | Memory (RAM) |
|--|----------------------------------|--------------------|
| Windows® 7 (64 bit) Professional with Service Pack 1 | 2.83 GHz Intel Pentium Quad Core | 16 GB ¹ |
| Windows® 10 (64 bit) Professional | 2.83 GHz Intel Pentium Quad Core | 16 GB ¹ |

¹ Systems with 8 GB can be used, however it is recommended you limit your processing to <200 samples per analyses.

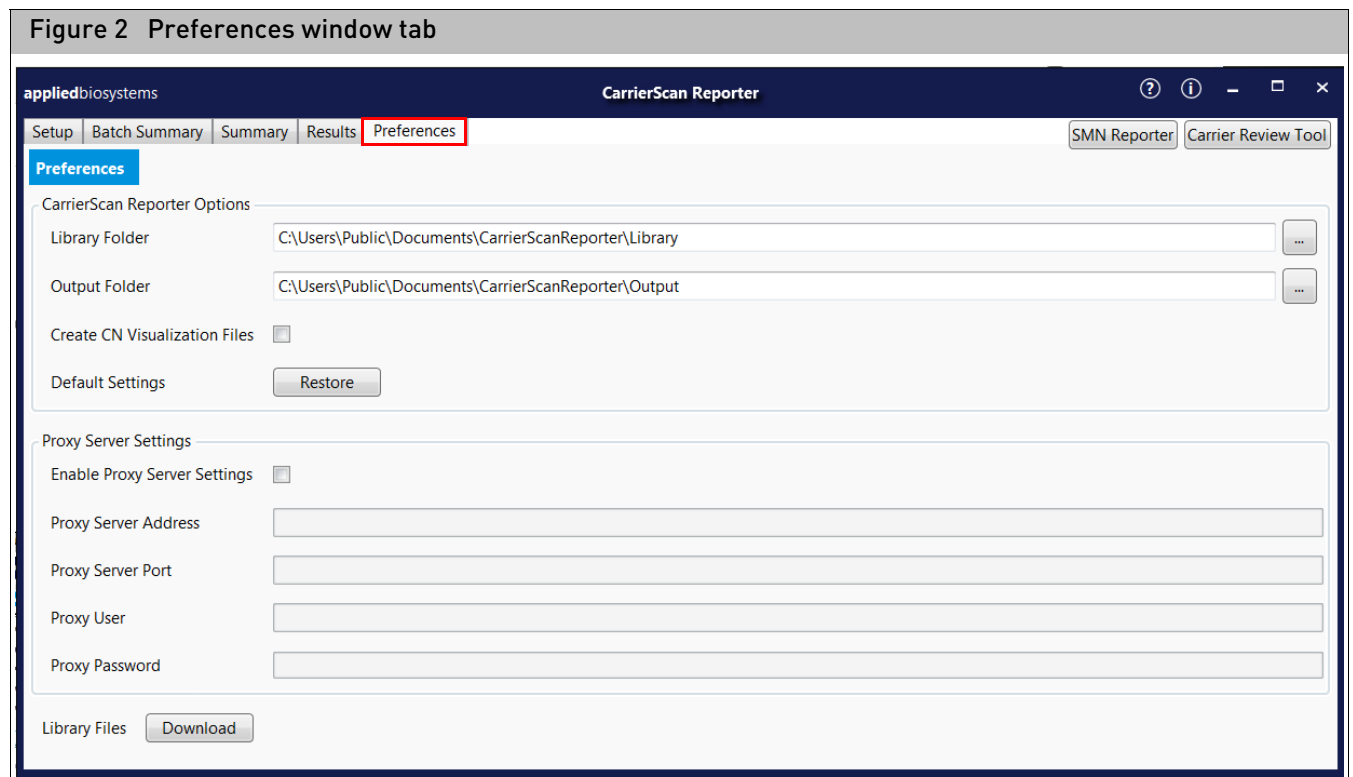
Installation

1. Go to: thermofisher.com
2. Locate and download the CarrierScan Reporter zipped software package.
3. Unzip the file as you normally would, then double-click on **CarrierScanReporterSetup.exe**.
4. Follow the installer's directions.
5. After the installation is complete, click **Start** → **All Programs** → **Thermo Fisher Scientific** → **CarrierScan Reporter**
The application opens.

Start up

The library and output folders are preset to default locations. Before using this software, review or change these folder paths, install required library files, and (if needed) set up access to your site's Proxy Server.

1. Click on the **Preferences** tab. (Figure 2)



Locating the library folder

The Library folder's default location is: C:\Users\Public\Documents\CarrierScanReporter\Library. If you want to change this default location, continue to Step 1.

1. Click the Library Folder field's **Browse** button, navigate to the location/folder you want, then click **Select Folder**.

Your assigned library folder path is now displayed.

Note: To better organize your data within your library folder, create a sub-folder for each supported array type.

Locating the output folder

This folder stores data your analysis generated. The Output folder's default location is: C:\Users\Public\Documents\CarrierScanReporter\Output. If you want to change this default location, continue to Step 1.

1. Click the Output Folder field's **Browse** button, navigate to the location/folder you want, then click **Select Folder**.

Your assigned output folder path is now displayed.

(Optional) Creating copy number visualization files

If you want to generate a *.cn file that contains log2 ratio data from all files in your analysis, click the **Create CN Visualization Files** check box. After the *.cn file is generated, it is saved to your assigned output folder.

This file can be viewed using an Integrative Genomics Viewer (IGV).

To access this viewer, go to: <http://software.broadinstitute.org/software/igv/>

Setting up custom proxy settings

Follow the steps below if your system has to pass through a Proxy Server before it can access the NetAffx server.

1. Click the **Enable Proxy Settings** check box.

Note: The proxy user ID and password is NOT the same ID and password used to connect to the NetAffx server.

2. Enter the **Address, Port, User, and Password**. Contact your IT department, if you do not know the proxy settings.

Downloading library files

1. Click the **Download** button.

The NetAffx User Information window appears.

2. Enter your NetAffx account email and password, then click **OK** or go to netaffx.com and click **Register** to sign up.

Note: If you are unable to connect to NetAffx, make sure you have an active Internet connection, and/or correct Proxy Server settings.

The NetAffx Update window appears.

3. Click the corresponding check box(es) of the library file(s) you want to download, then click **OK**.

An Installing Updates progress bar appears.

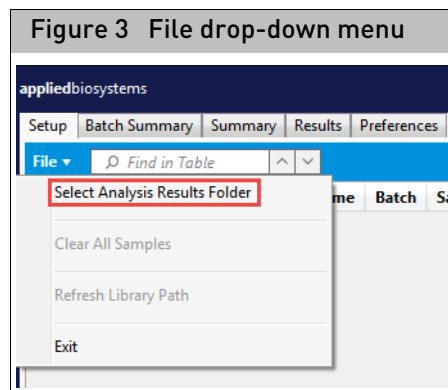
4. Click **OK** to close the NetAffx Update window.

Note: If creating a custom annotation panel file, see *Note* in the full TXT report.

IMPORTANT! Before you can analyze data using CarrierScan Reporter, your CEL files MUST first be processed through the Axiom Analysis Suite 3.1 or higher software application. For instructions on how to do this, refer to the Axiom Analysis Suite User Guide (P/N 703307).

Importing data

1. Click **File** → **Select Analysis Results Folder**. (Figure 3)



A window opens.

2. Navigate to the Axiom Analysis Suite output folder.
3. Open the Analysis Results Folder that has the samples you want to analyze.
4. Click **Select Folder**.

Note: Multiple Analysis Results Folders can be added into the same analysis. To do this, repeat steps 1-4.

IMPORTANT! Analysis Results Folder names must be unique and subsequently loaded results folders must be for the same array type and analysis settings as the first folder.

The sample files from your selected Analysis Results Folder(s) load into the Setup window tab. (Figure 4)

Note: The Setup table (Figure 4) may take longer to populate if you are loading a large file and/or checked the **Create CN Visualization Files** check box (in Preferences).

Figure 4 Loaded data in the Setup window tab

| Panel | Analysis Name | Batch | Sample Name | M/F | DQC | QC Call Rate | MAPD |
|-------|---------------|-----------------------------------|--|-----|-------|--------------|-------|
| | | mPCR102_P2_nostep2.cn.v3_20181009 | CD00021_CscanTraining_FX_ON_P2_Wk1_C07 | M | 0.993 | 100.0 | 0.127 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA00006_CscanTraining_FX_ON_P2_Wk1_H08 | F | 0.997 | 100.0 | 0.139 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA00937_CscanTraining_FX_ON_P2_Wk1_A09 | M | 0.990 | 100.0 | 0.143 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA01607_CscanTraining_FX_ON_P2_Wk1_F04 | M | 0.994 | 100.0 | 0.130 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA02533_CscanTraining_FX_ON_P2_Wk1_D05 | F | 0.993 | 100.0 | 0.143 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA05046_CscanTraining_FX_ON_P2_Wk1_C05 | F | 0.995 | 100.0 | 0.136 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA05117_CscanTraining_FX_ON_P2_Wk1_H06 | F | 0.993 | 99.9 | 0.139 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA05159_CscanTraining_FX_ON_P2_Wk1_G06 | F | 0.991 | 100.0 | 0.130 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA05816_CscanTraining_FX_ON_P2_Wk1_B11 | F | 0.995 | 100.0 | 0.157 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA08684_CscanTraining_FX_ON_P2_Wk1_C11 | F | 0.996 | 100.0 | 0.129 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA11468_CscanTraining_FX_ON_P2_Wk1_D09 | F | 0.997 | 100.0 | 0.143 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA11472_CscanTraining_FX_ON_P2_Wk1_A02 | M | 0.996 | 100.0 | 0.131 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA11723_CscanTraining_FX_ON_P2_Wk1_B02 | F | 0.993 | 100.0 | 0.146 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA11761_CscanTraining_FX_ON_P2_Wk1_H01 | M | 0.995 | 100.0 | 0.139 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA12960_CscanTraining_FX_ON_P2_Wk1_E03 | M | 0.995 | 100.0 | 0.138 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA13591_CscanTraining_FX_ON_P2_Wk1_E02 | F | 0.995 | 100.0 | 0.135 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA14108_CscanTraining_FX_ON_P2_Wk1_F09 | F | 0.995 | 100.0 | 0.126 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA16266_CscanTraining_FX_ON_P2_Wk1_F07 | M | 0.993 | 99.8 | 0.140 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA17912_CscanTraining_FX_ON_P2_Wk1_E08 | F | 0.996 | 100.0 | 0.131 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA18445_CscanTraining_FX_ON_P2_Wk1_C08 | F | 0.992 | 100.0 | 0.142 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA18929_CscanTraining_FX_ON_P2_Wk1_A01 | M | 0.996 | 99.9 | 0.170 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA20945_CscanTraining_FX_ON_P2_Wk1_E11 | F | 0.993 | 100.0 | 0.134 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA21081_CscanTraining_FX_ON_P2_Wk1_F06 | M | 0.998 | 99.9 | 0.147 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA23687_CscanTraining_FX_ON_P2_Wk1_A06 | F | 0.995 | 100.0 | 0.144 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA20942_CscanTraining_FX_ON_P2_Wk1_D11 | M | 0.995 | 99.9 | 0.131 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA21551_CscanTraining_FX_ON_P2_Wk1_D01 | F | 0.997 | 100.0 | 0.143 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA23087_CscanTraining_FX_ON_P2_Wk1_A07 | F | 0.990 | 100.0 | 0.132 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA23437_CscanTraining_FX_ON_P2_Wk1_A12 | F | 0.993 | 99.9 | 0.140 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | CD00023_CscanTraining_FX_ON_P2_Wk1_D07 | F | 0.994 | 100.0 | 0.132 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA00449_CscanTraining_FX_ON_P2_Wk1_B06 | F | 0.996 | 100.0 | 0.142 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA00879_CscanTraining_FX_ON_P2_Wk1_G11 | F | 0.996 | 100.0 | 0.136 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA02795_CscanTraining_FX_ON_P2_Wk1_E10 | M | 0.993 | 100.0 | 0.130 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA03252_CscanTraining_FX_ON_P2_Wk1_F05 | F | 0.998 | 100.0 | 0.142 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA04408_CscanTraining_FX_ON_P2_Wk1_C01 | M | 0.998 | 99.8 | 0.138 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA10799_CscanTraining_FX_ON_P2_Wk1_E06 | M | 0.995 | 100.0 | 0.136 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA11275_CscanTraining_FX_ON_P2_Wk1_D03 | M | 0.996 | 99.9 | 0.140 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA17436_CscanTraining_FX_ON_P2_Wk1_B10 | M | 0.995 | 99.9 | 0.144 |
| | | mPCR102_P2_nostep2.cn.v3_20181009 | NA04258_CscanTraining_FX_ON_P2_Wk1_A11 | M | 0.994 | 99.9 | 0.145 |

You can sort the data (in ascending or descending order) in each column by clicking on its header. See the table below for column definitions.

Note: Once an analysis has been run and samples are loaded into the table, a QC file is written as **SampleQC.txt**, where all QC values shown on the Setup table are recorded for all samples analyzed.

| Column | Description |
|---------------|---|
| Panel | Displays the name of the panel used in the analysis. |
| Analysis Name | Analysis Name provided in the Analysis Mode or the Mapping File. |
| Batch | Name assigned to the collection of data in the selected analysis results folder. |
| Sample Name | The assigned sample name (that was designated during its sample registration process in Command Console). |
| M/F | M = Male, F = Female, U = Unknown |
| DQC (Dish QC) | DQC is based on intensities of probe sequences for non-polymorphic genome locations (i.e., sites that do not vary in sequence from one individual to the next). When subject to the two-color Axiom assay, probes expected to ligate an A or T base (referred to as AT non-polymorphic probes) produce specific signal in the AT channel and background signal in the GC channel. The converse is true for probes expected to ligate a G or C base (referred to as GC non-polymorphic probes). DQC is a measure of the resolution of the distributions of "contrast" values, where: Distributions of contrast values are computed separately for the AT non-polymorphic probes (which should produce positive contrast values) and GC non-polymorphic probes (which should produce negative contrast values). If sample quality is high, then signal will be high in the expected channel and low in background channel, and the two contrast distributions will be well-resolved. A DQC value of zero indicates no resolution between the distributions of AT and GC probe contrast values, and the value of 1 indicates perfect resolution. |
| QC Call Rate | Percentage of SNPs assigned a genotype using a subset of probe sets (usually 20,000) that are autosomal. |
| MAPD | Median of the Absolute values of all Pairwise Differences (MAPD) is a global measure of the variation of all microarray probes across the genome. It represents the median of the distribution of changes in Log2 Ratio between adjacent probes. Since it measures differences between adjacent probes, it is a measure of short-range noise in the microarray data. |
| MAPDc | MAPD calculation post plate correction. |
| WavinessSD | A global measure of variation of microarray probes that is insensitive to short-range variation and focuses on long-range variation. |
| WavinessSDc | WavinessSD calculation post plate correction. |

Selecting an analysis mode

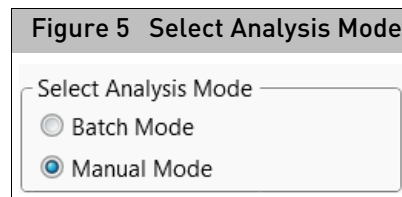
You can setup your analysis in two ways:

- **Manual Mode** supports individual sample analysis set up, paired sample analysis, and modifications. In this mode, samples can be added one by one or as pairs.
- **Batch Mode** supports quick, pre-determined setup analysis through a structured mapping file.

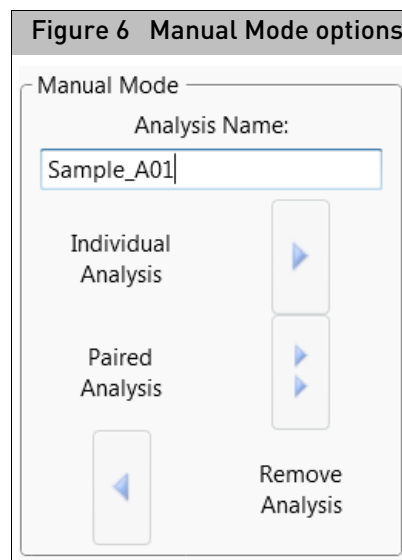
Either mode enables you to select and run one analysis at a time.

Manual mode

1. Click the **Manual Mode** radio button. (Figure 5)



The Manual Mode Options window appears. (Figure 6)

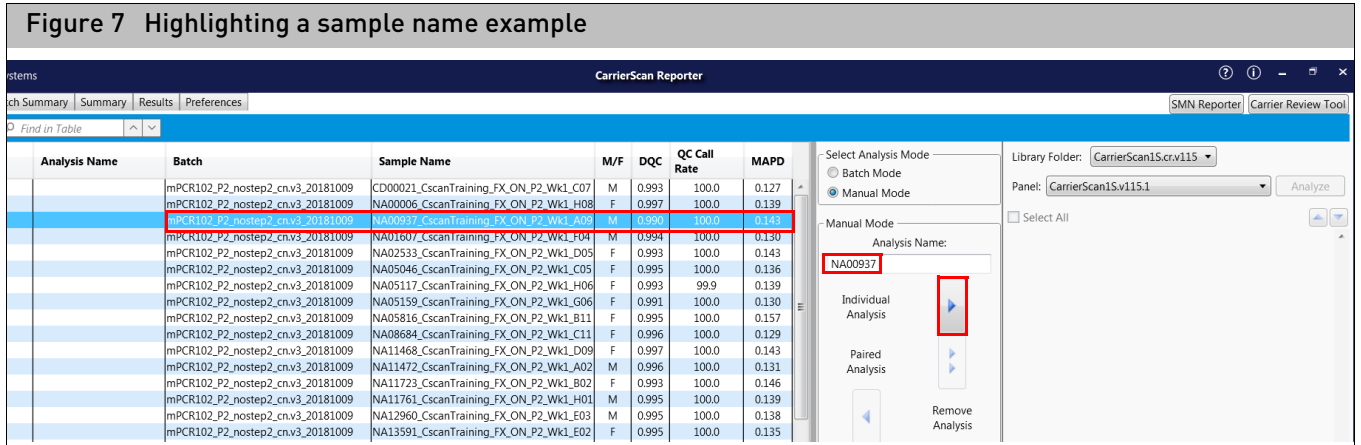


Individual analysis (one sample)

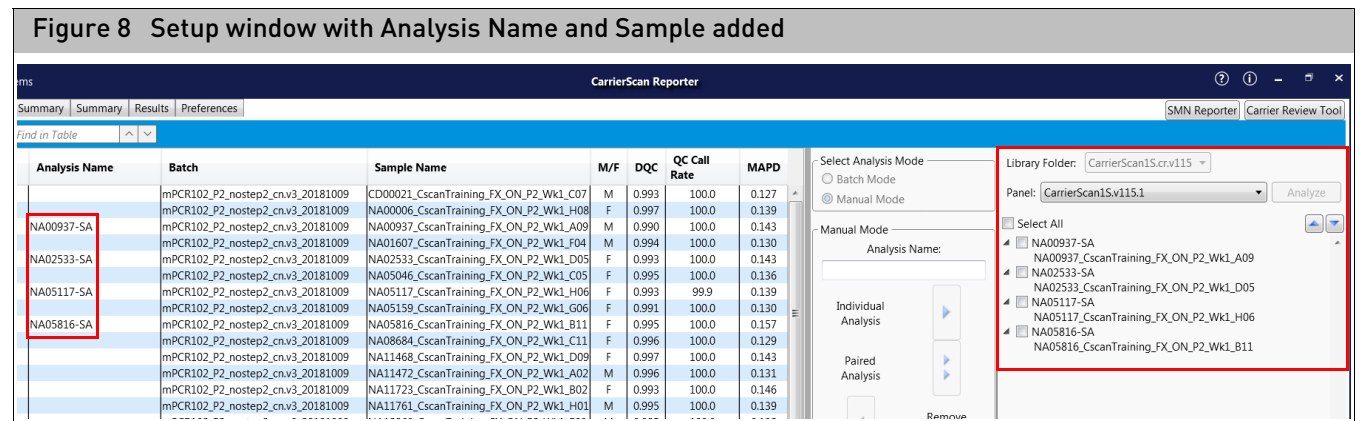
1. Highlight a sample name in the table, as shown in Figure 7.
2. Enter an Analysis Name. If no Analysis Name is provided, the sample name will be used as the Analysis Name.

Note: "SA" is appended to the analysis name to indicate the analysis is a Single Analysis.

- Click the newly enabled **Individual Analysis** arrow button (Figure 7) to add the sample to the Analysis Window.



The Analysis Name and its associated sample are now added to the far right-hand side of the Analysis Window. The Analysis Name is also added to the Table's **Analysis Name** column, as shown in Figure 8.



Adding multiple samples to an individual analysis simultaneously

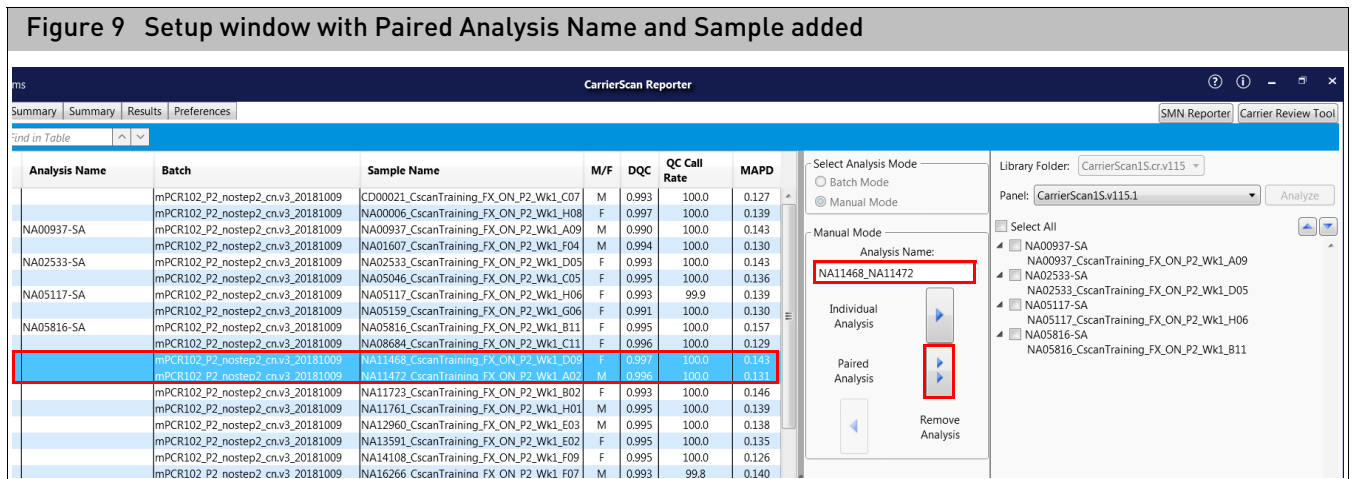
- Highlight the desired sample names in the table.
- Enter an Analysis Name. If no Analysis Name is provided, the sample name will be used as the Analysis Name.
- Click the newly enabled Individual Analysis arrow to add the sample to the Analysis Window.

The Analysis Name and its associated sample are added to the far right-hand side of the Analysis Window and are now ready for analysis. The Analysis Name is also added to the Table's **Analysis Name** column, as shown in Figure 8.

Paired analysis (two samples)

1. Click on your first sample, then press the Ctrl key and click on your second sample.
Both samples are highlighted, as shown in Figure 9.
2. Enter an Analysis Name. If an Analysis Name is not provided, the name of the first sample highlighted for the paired analysis will be used as the Analysis Name.
Note: PA is appended to the analysis name to indicate the analysis is a Paired Analysis.
3. Click the newly enabled Paired Analysis arrow to add the sample to the Analysis Window, as shown in Figure 9.

Figure 9 Setup window with Paired Analysis Name and Sample added



Removing an analysis name from the setup

1. Check the check box of the Analysis Name you want to remove.
2. Click on the newly enabled **Remove Analysis** button.

Running an analysis

1. Select the appropriate panel from the Panel drop-down.
2. Check the Analysis Name(s) to be analyzed with that panel or click the **Check All** box to include all analysis names in the analysis.
3. Once all analyses have been added, click the **Analyze** button.

IMPORTANT! You can only add one sample per analysis. If you have added a sample that already exists in the analysis you want to run, a window appears detailing why your sample could not be added. Acknowledge the message, then click **OK**.

The analysis process begins and a Please Wait window appears.

After the analysis is complete, the remaining columns in the table are populated, as shown in Figure 10.

Figure 10 Populated Setup window

The screenshot displays the CarrierScan Reporter application window. The main area contains a table with the following columns: Analysis Name, Batch, Sample Name, M/F, DQC, QC Call Rate, and MAPD. The table is populated with numerous rows of data, including sample IDs like CD00021, NA00006, and NA00937, and their corresponding QC metrics.

On the right side of the window, there is a control panel with the following sections:

- Select Analysis Mode:** Radio buttons for Batch Mode and Manual Mode.
- Manual Mode:** A text field for Analysis Name and three buttons: Individual Analysis, Paired Analysis, and Remove Analysis.
- Library Folder:** A dropdown menu set to CarrierScan1S.cr.v115.1.
- Panel:** A dropdown menu set to CarrierScan1S.v115.1 and an Analyze button.
- Select All:** A checkbox.
- Tree View:** A hierarchical list of analysis items, including folders for NA00937-SA, NA02533-SA, NA05117-SA, NA05816-SA, and NA11468_NA11472-PA, with sub-items for various sample names.

Batch mode

IMPORTANT! Before using the Batch Mode feature, you must first create a tab-delimited TXT file (Figure 11) to define your individual or paired analysis.

Required batch mode file guidelines

Your Batch Mode Mapping file must contain these six header columns (Figure 11) in the order shown:

[A] Panel - The name of the panel for the analysis.

[B] Analysis Name - The name given to the analysis of that sample(s). **Note:** For a paired analysis, use the same analysis name for both samples in the pair.

[C] Batch - This name must match the batch name provided when loading the samples.

[D] Sample Name - The filename of the samples to be analyzed.

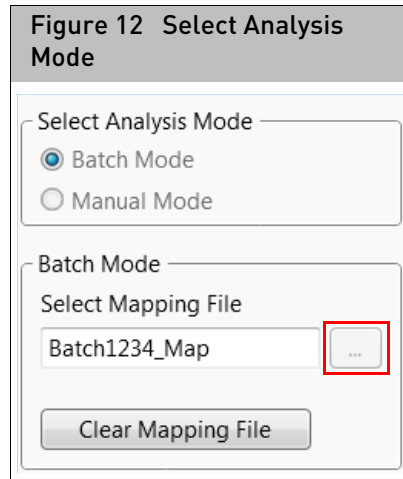
[E] Ethnicity - The assignment of ethnicity only applies to paired analyses. If a valid ethnicity (present on the chosen panel) is assigned to one sample, then the paired sample must also have a valid ethnicity assigned. Ethnicities are used in the pair residual chance of occurrence calculations provided in the full and positive outputs. Ethnicity assignments are not required. **Note:** Column [E] is optional. If you choose not to include it, use your space bar to insert a single blank space in each of its cells to ensure a properly formatted TSV file.

[F] Notes - Use this column to enter any comments you want regarding the samples or analysis. Cells in this column can be left blank, as comments are not required. **Note:** Column [F] is optional. If you choose not to include it, use your space bar to insert a single blank space in each of its cells to ensure a properly formatted TSV file.

Figure 11 Required and optional tab-delimited TXT file example

| | A | B | C | D | E | F |
|----|---------------------|-----------------|---------------------|-------------|--------------------|-----------|
| 1 | Panel | Analysis Name | Batch | Sample Name | Ethnicity | Notes |
| 2 | CarrierScan.r1 | CD00008 | TrainingPlate_r1_EX | CD00008_C02 | | expedite |
| 3 | CarrierScan.r1.EX | CD00021_CD00023 | TrainingPlate_r1_EX | CD00021_C07 | General Population | paired |
| 4 | CarrierScan.r1.EX | CD00021_CD00023 | TrainingPlate_r1_EX | CD00023_D07 | Asian American | paired |
| 5 | CarrierScan.r1.CFTR | Run180103_H08 | TrainingPlate_r1_EX | NA00006_H08 | | CFTR only |
| 6 | CarrierScan.r1 | Sample_NA00059 | TrainingPlate_r1_EX | NA00059_A04 | | |
| 7 | CarrierScan.r1.ACOG | NA00244_G09 | TrainingPlate_r1_EX | NA00244_G09 | | ACOG only |
| 8 | CarrierScan.r1 | NA00449_B06 | TrainingPlate_r1_EX | NA00449_B06 | | rerun |
| 9 | CarrierScan.r1.CFTR | Pair-E09_E04 | TrainingPlate_r1_EX | NA00649_E09 | Caucasian | |
| 10 | CarrierScan.r1.CFTR | Pair-E09_E04 | TrainingPlate_r1_EX | NA00852_E04 | Caucasian | |

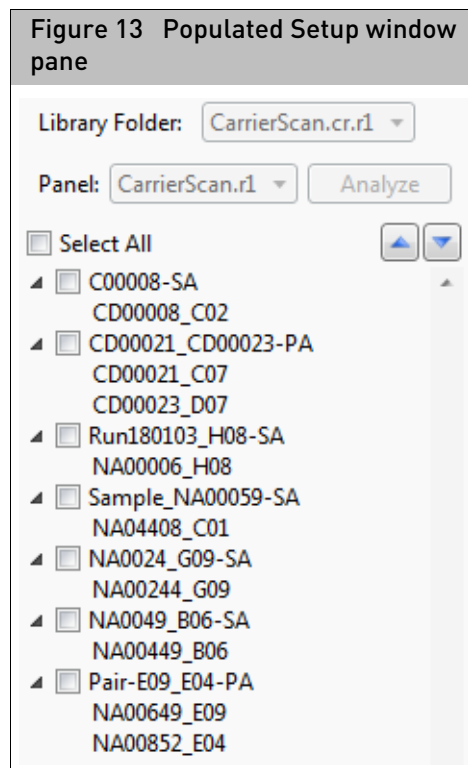
1. After creating and saving your tab-delimited Batch Mode TXT file, click the **Batch Mode** radio button. (Figure 12)



If you want to remove samples from the Analysis window, click the **Clear Mapping File** button.

2. Click **Select Mapping File** Browse button.
An Explorer window appears.
3. Navigate to the Batch Mode file, then click **OK**.

The Analysis Setup window populates with individual and paired analysis (based on the information in your Batch Mode file), as shown in Figure 13.



- Use the check box to check each Analysis Name(s) to be analyzed or click the **Check All** check box to include all the names shown in your analysis.
- Click the **Analyze** button.

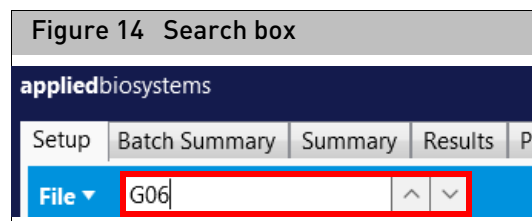
The analysis process begins and a Please Wait window appears. After the analysis is complete, the remaining columns in the table are populated.

Searching table content

Use this feature to search for batch, gender, and sample names within the table.

Note: Every column, except for the Analysis Completed column is searched.

- From the Analysis Setup window tab's **Search** box (Figure 14), type the name or portion of the name you want to find.



- Click the up or down arrow (or press the Enter key).

If a match is found, it is highlighted in the table, as shown in Figure 15.

Figure 15 Found (highlighted) entry

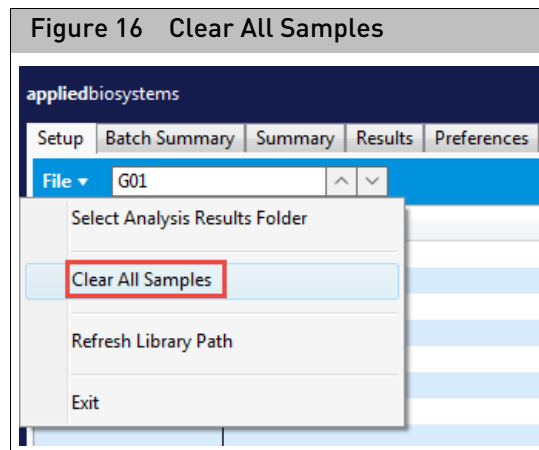
| appliedbiosystems | | | | | | | | |
|----------------------|--------------------|-----------------------------------|--|-------------|-------|--------------|-------|--|
| CarrierScan Reporter | | | | | | | | |
| Setup | Batch Summary | Summary | Results | Preferences | | | | |
| File | G06 | | | | | | | |
| Panel | Analysis Name | Batch | Sample Name | M/F | DQC | QC Call Rate | MAPD | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | CD00021_CscanTraining_FX_ON_P2_Wk1_C07 | M | 0.993 | 100.0 | 0.127 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA00006_CscanTraining_FX_ON_P2_Wk1_H08 | F | 0.997 | 100.0 | 0.139 | |
| | NA00937-SA | mPCR102_P2_nostep2_cn.v3_20181009 | NA00937_CscanTraining_FX_ON_P2_Wk1_A09 | M | 0.990 | 100.0 | 0.143 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA01607_CscanTraining_FX_ON_P2_Wk1_F04 | M | 0.994 | 100.0 | 0.130 | |
| | NA02533-SA | mPCR102_P2_nostep2_cn.v3_20181009 | NA02533_CscanTraining_FX_ON_P2_Wk1_D05 | F | 0.993 | 100.0 | 0.143 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA05046_CscanTraining_FX_ON_P2_Wk1_C05 | F | 0.995 | 100.0 | 0.136 | |
| | NA05117-SA | mPCR102_P2_nostep2_cn.v3_20181009 | NA05117_CscanTraining_FX_ON_P2_Wk1_H06 | F | 0.993 | 99.9 | 0.139 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA05159_CscanTraining_FX_ON_P2_Wk1_G06 | F | 0.991 | 100.0 | 0.130 | |
| | NA05816-SA | mPCR102_P2_nostep2_cn.v3_20181009 | NA05816_CscanTraining_FX_ON_P2_Wk1_B11 | F | 0.995 | 100.0 | 0.157 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA08684_CscanTraining_FX_ON_P2_Wk1_C11 | F | 0.996 | 100.0 | 0.129 | |
| | NA11468_NA11472-PA | mPCR102_P2_nostep2_cn.v3_20181009 | NA11468_CscanTraining_FX_ON_P2_Wk1_D09 | F | 0.997 | 100.0 | 0.143 | |
| | NA11468_NA11472-PA | mPCR102_P2_nostep2_cn.v3_20181009 | NA11472_CscanTraining_FX_ON_P2_Wk1_A02 | M | 0.996 | 100.0 | 0.131 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA11723_CscanTraining_FX_ON_P2_Wk1_B02 | F | 0.993 | 100.0 | 0.146 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA11761_CscanTraining_FX_ON_P2_Wk1_H01 | M | 0.995 | 100.0 | 0.139 | |
| | | mPCR102_P2_nostep2_cn.v3_20181009 | NA12960_CscanTraining_FX_ON_P2_Wk1_E03 | M | 0.995 | 100.0 | 0.138 | |

- Click the up or down arrow (or press the Enter key) to go to the next/previous matching entry.

Removing all samples from the table

1. Click **File** → **Clear All Samples**. (Figure 16)
All Samples are removed from the table.

IMPORTANT! Any unanalyzed analyses (shown in the right panel) are also removed.

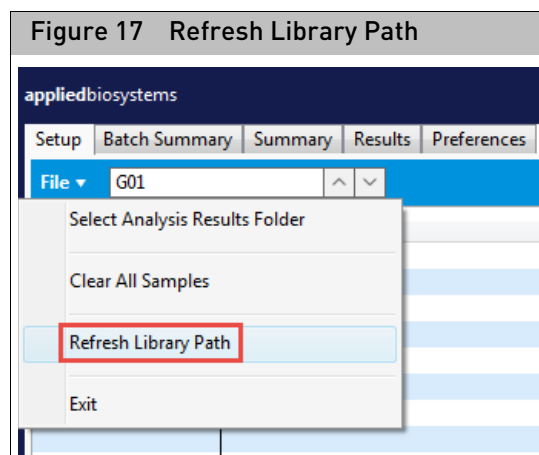


Updating the library folder

To update the contents of your library folder:

1. Click **File** → **Refresh Library Path**. (Figure 17)

The Library folder is now updated to include your changes (since the application was last opened) and any available library file updates.



3

Summary tables

Batch summary table

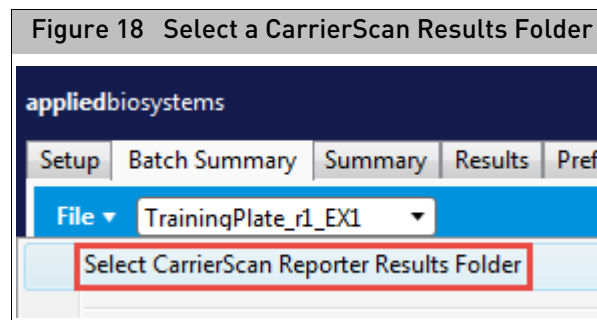
After setting up and successfully running analyses for an Axiom Analysis Suite Results Folder as described in [Chapter 2, "Setup"](#), click the **Batch Summary** tab to visually inspect variants across all samples analyzed within an Axiom Analysis Suite Results Folder (also referred to as the Batch).

Note: If multiple panels are used with a batch, each panel will generate its own BatchSummaryResults file.

The Batch Summary table contains information for all samples analyzed in a batch. Viewing results for a 96 sample plate is supported in the software, however we advise to view results >96 samples in other software (e.g. Excel) to minimize slowing of software performance.

Selecting previously generated results

1. From the **Batch Summary** window tab, click **File** → **Select CarrierScan Reporter Results Folder**. ([Figure 18](#))



A **Select results folder** window appears.

2. Click to highlight the analysis folder you want to view in detail, then click **Select Folder**.

Note: Past results can only be viewed if the original files (inside the folder you select) have not been modified.

A detailed analysis of your selected folder appears, as shown in [Figure 19](#).

Figure 19 Batch Summary table

| # | Gene:Variant | 5509544367216051220568_(CarrierScanIS)_A01 (4) | 5509544367216051220568_(CarrierScanIS)_A02 (3) | 5509544367216051220568_(CarrierScanIS)_A03 (3) | 5509544367216051220568_(CarrierScanIS)_A04 (3) | 5509544367216051220568_(CarrierScanIS)_A05 (2) | 5509544367216051220568_(CarrierScanIS)_A06 (4) | 5509544367216051220568_(CarrierScanIS)_A07 (4) | 5509544367216051220568_(CarrierScanIS)_A08 (3) | 5509544367216051220568_(CarrierScanIS)_A09 (3) | 5509544367216051220568_(CarrierScanIS)_A10 (4) | 5509544367216051220568_(CarrierScanIS)_A11 (4) | 5509544367216051220568_(CarrierScanIS)_A12 (2) | 5509544367216051220568_(CarrierScanIS)_B01 (3) | 5509544367216051220568_(CarrierScanIS)_B02 (3) | 5509544367216051220568_(CarrierScanIS)_B03 (2) | 5509544367216051220568_(CarrierScanIS)_B04 (3) | 5509544367216051220568_(CarrierScanIS)_B05 (3) |
|---|---|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|
| 2 | CFTR:c.1647T>G pathogenic; drug response | N | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | SERPINA1:c.1096G>A pathogenic; other | N | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.1865G>A pathogenic | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.3718-3T>G pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | BCHEC:c.293A>G pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.1364C>A pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | ATP7B:c.2333G>T pathogenic | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.3472C>T pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | HEXA:c.533G>A pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | CYP21A2:c.1174G>A likely pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | DHCR7:c.964-1G>C pathogenic/likely pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | CYP17A1:c.1084C>T pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | DNAH5:c.13486C>T pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | FAH:c.1062+5G>A pathogenic | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |

Click on a column header to sort the data (in ascending or descending order). See the table below for column definitions.

IMPORTANT! As additional samples are run, you must refresh the Batch Summary table to reveal the newly analyzed samples. To do this, click **File** → **Select CarrierScan Results Folder**.

| Column | Description |
|----------------|---|
| # | Number of samples with Carrier (C), Affected (A), or CNgain statuses. |
| Gene Variant | Identity of the Variant as the combination of the Gene and Variant Status as found on the panel in their respective columns. Note: Clicking on the Gene:Variant auto-launches a ClinVar search result for the Variant (if it exists). For exonic regions, you must perform the ClinVar search manually. |
| Sample Columns | Sample Names with the number of variants with Carrier (C), Affected (A), Unaffected Normal (N), or CNgain statuses. Note: A sample status with an asterisk [*] denotes multiple sample status. |

Selecting a sample column will auto-sort the column by its sample status (Figure 20) in the following order:

- Affected (A)
- Carrier (C)
- CN gain
- NoCall/Unknown (NC)
- NRP
- Normal (N)

Figure 20 Sample sorting

| | | NA11723 | NA11761 | NA11859 | NA12585 | NA12785 | NA12794 | NA12960 | NA12961 | NA13205 | NA13423 | NA13591 | NA14108 | NA16028 | NA16193 | NA16266 | NA17431 | NA17436 | NA17821 | NA17912 | NA18397 | NA18445 | NA18668 | NA18800 | NA18929 | NA20270 | NA20732 | NA20744 | NA20847 | NA20915 | NA20925 |
|----|---|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|---------|
| 54 | MTHFR:c.665C>T[conflicting] | A | N | A | C | N | C | C | N | C | C | C | C | C | C | N | C | C | A | C | N | N | N | N | N | N | A | N | N | C | C |
| 4 | CFTR:c.1210-7_1210-6delTT[Unknown Significance] | A | N | N | N | N | N | N | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.3846G>A[unknown] | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | C | N | N | N | N | N |
| 47 | ASPA:c.693C>T[benign] | C | N | N | N | N | C | C | N | C | C | N | N | N | C | N | N | N | C | A | C | C | C | C | N | C | N | N | A | C | N |
| 1 | HEXA:c.1421+1G>C[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HEXA:c.1274_1277dupTATC[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | HEXA:c.1073+1G>A[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 0 | HEXA:c.-2564[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HBB:c.92+6T>C[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 0 | HBB:c.79G>T[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HBB:c.79G>A[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HBB:c.316-149[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HBB:c.316-106C>G[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 1 | HRR:r.20rdeA[unknown] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |

Displaying only pathogenic variants

1. Click the **Show only pathogenic variants** check box.

Only variants with pathogenic statuses from the panel appear.

Note: Sample tallies within the column headers are recalculated based on what is displayed, however the # column does NOT recalculate to show which samples have Carrier (C), Affected (A) or CNgain statuses for that variant, as shown in Figure 21.

Figure 21 Show only pathogenic variants

| # | Gene:Variant | 5509544367216051220568_(CarrierScan15)_A01 (4) | 5509544367216051220568_(CarrierScan15)_A02 (3) | 5509544367216051220568_(CarrierScan15)_A03 (3) | 5509544367216051220568_(CarrierScan15)_A04 (3) | 5509544367216051220568_(CarrierScan15)_A05 (2) | 5509544367216051220568_(CarrierScan15)_A06 (4) | 5509544367216051220568_(CarrierScan15)_A07 (4) | 5509544367216051220568_(CarrierScan15)_A08 (3) | 5509544367216051220568_(CarrierScan15)_A09 (3) | 5509544367216051220568_(CarrierScan15)_A10 (4) | 5509544367216051220568_(CarrierScan15)_A11 (4) | 5509544367216051220568_(CarrierScan15)_A12 (2) | 5509544367216051220568_(CarrierScan15)_B01 (3) | 5509544367216051220568_(CarrierScan15)_B02 (3) | 5509544367216051220568_(CarrierScan15)_B03 (2) | 5509544367216051220568_(CarrierScan15)_B04 (3) | 5509544367216051220568_(CarrierScan15)_B05 (3) | 5509544367216051220568_(CarrierScan15)_B06 (3) | 5509544367216051220568_(CarrierScan15)_B07 (4) | 5509544367216051220568_(CarrierScan15)_B08 (4) |
|---|---|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|--|
| 2 | CFTR:c.1647T>G[pathogenic; drug response] | N | N | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | SERPINA1:c.1096G>A[pathogenic; other] | N | N | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.1865G>A[pathogenic] | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.3718-3T>G[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | BCHEC:c.293A>G[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | CFTR:c.1364C>A[pathogenic] | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |
| 2 | ATP7B:c.2333G>T[pathogenic] | N | N | N | C | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N | N |

Summary table

After setting up and successfully running an analysis, as described in [Chapter 2](#), "Setup" on page 9, click the **Summary** window tab to view your results. (Figure 22)

The Summary table provides a summary of all analyses run in a session. Each session's analyses are contained within a time-stamped folder. This folder includes the first session's analysis and any subsequent analyses thereafter - until the software is closed.

An **AnalysisSummary.txt** file is auto-generated at the start of each session. This file captures your session's table data, then continues to auto-save any subsequent analyses until you end your session. The file resides in the Output folder you assigned earlier and can be used for reference purposes.

Figure 22 Summary window tab

| Analysis Name | Sample Name | Batch Name | Sample RC3 | Coupled Sample Name | Coupled Sample Batch Name | Coupled Sample RC3 | Variants Detected | Variants Tested | Associated Genes | Associated Phenotypes | Common Marked Genes | Variants To Review | Panel Analyzed |
|---|--|----------------------------------|------------|--|----------------------------------|--------------------|-------------------|-----------------|------------------|-----------------------|---------------------|--------------------|------------------|
| CD00023_CscanTraining_FX_ON_P2_Wk1_D07-SA | CD00023_CscanTraining_FX_ON_P2_Wk1_D07 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 13 | 5667 | 10 | 7 | | 4 | Draft.20190403.r |
| NA00449_CscanTraining_FX_ON_P2_Wk1_B06-SA | NA00449_CscanTraining_FX_ON_P2_Wk1_B06 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 13 | 5666 | 11 | 8 | | 5 | Draft.20190403.r |
| NA00879_CscanTraining_FX_ON_P2_Wk1_G11-SA | NA00879_CscanTraining_FX_ON_P2_Wk1_G11 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 12 | 5666 | 11 | 8 | | 5 | Draft.20190403.r |
| NA02795_CscanTraining_FX_ON_P2_Wk1_E10-SA | NA02795_CscanTraining_FX_ON_P2_Wk1_E10 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 14 | 5659 | 13 | 9 | | 6 | Draft.20190403.r |
| NA03252_CscanTraining_FX_ON_P2_Wk1_F05-SA | NA03252_CscanTraining_FX_ON_P2_Wk1_F05 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 12 | 5667 | 12 | 9 | | 5 | Draft.20190403.r |
| NA04258_CscanTraining_FX_ON_P2_Wk1_A11-SA | NA04258_CscanTraining_FX_ON_P2_Wk1_A11 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 12 | 5667 | 11 | 7 | | 6 | Draft.20190403.r |
| NA04395_CscanTraining_FX_ON_P2_Wk1_H10-PA | NA04395_CscanTraining_FX_ON_P2_Wk1_H10 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | NA05258_CscanTraining_FX_ON_P2_Wk1_F08 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | 24 | 11331 | 17 | 11 | 1 | 7 | Draft.20190403.r |
| NA04408_CscanTraining_FX_ON_P2_Wk1_C01-SA | NA04408_CscanTraining_FX_ON_P2_Wk1_C01 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | | | | 14 | 5667 | 13 | 8 | | 7 | Draft.20190403.r |
| NA07441_CscanTraining_FX_ON_P2_Wk1_G03-PA | NA07441_CscanTraining_FX_ON_P2_Wk1_G03 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | NA11067_CscanTraining_FX_ON_P2_Wk1_G03 | mPCR102_P2_nostep2_cnv3_20181009 | Pass | 24 | 11333 | 13 | 8 | 2 | 6 | Draft.20190403.r |

Click on a column header to sort the data (in ascending or descending order). See the table below for column definitions.

| Column | Description |
|---------------------------|---|
| Analysis Name | The name provided for the analysis. SA denotes individual analysis. PA denotes paired analysis. |
| Sample Name | Displays the filename of the sample. |
| Batch Name | The name provided for the Batch that contains the sample. |
| Sample RC3 | Sample QC metric based on the number of rare variants determined to be non-normal. |
| Coupled Sample Name | Only for paired analysis. The other sample filename in the pair. |
| Coupled Sample Batch Name | Only for paired analysis. The Batch name containing the other filename in the pair. |
| Variants Detected | Number of non-normal genetic variations detected. |
| Variants Tested | Total genetic variations analyzed on the selected panel. |
| Associated Genes | Number of genes associated with the number of genetic variations detected. |
| Associated Phenotypes | Number of phenotypes associated with the number of genetic variations detected. |
| Panel Analyzed | Name of the Panel used in the analysis. |

| Column | Description |
|---------------------|--|
| Common Marked Genes | Number of common genes with 'pathogenic' genetic aberrations for paired samples. Pathogenic variants are defined by the Variant Status from the panel used for analysis. |
| Variants to Review | Number of genetic variants on the panel that require follow-up and review. Variants in this category were either Unknown or No Call. |
| Notes | <p>Notes that were provided on the Mapping File. Additional notes can be added.</p> <p><i>To do this:</i></p> <ol style="list-style-type: none"> 1. Click inside a Notes cell. A cursor appears. 2. Enter any additional notes. 3. Click outside the cell to add/save your additional note. <p>Note: Your amended notes are not saved unless you click File → Export and save the file using a different name.</p> |

Searching table content

1. From the Summary window tab's **Search** box, type the name or portion of the name you want to find.
2. Click the up or down arrow (or press the Enter key).
If a match is found, it is highlighted in the table.
3. Click the up or down arrow (or press the Enter key) to go to the next/previous matching entry.

Note: All columns are searched, except for columns containing date and/or timestamps.

Exporting the summary table

1. Click **File** → **Export Summary Results**.
An Explorer window appears.
2. Navigate to an easily accessible save location, name your export file, then click **Save**.

Your exported file is saved as a tab-delimited TXT file.

4

Results table

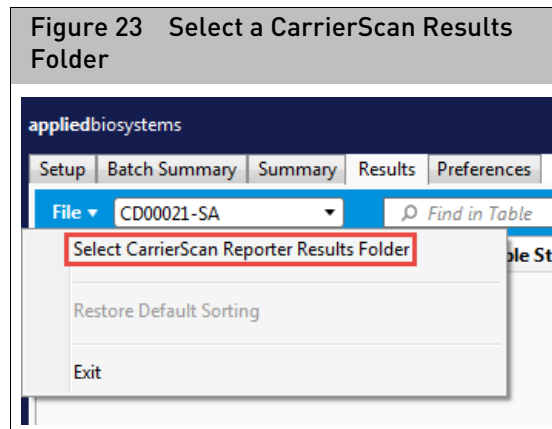
Using the results table

After reviewing your summary, as described in [Chapter 3, "Summary tables"](#), you can view more detailed results of your analysis by clicking on the **Results** window tab.

Note: If an analysis has been run in the current session, the default Results view lists your results by Analysis Name (sorted by lowest alphabetical name first).

Selecting a results folder

1. From the **Results** window tab, click **File** → **Select CarrierScan Reporter Results Folder**. ([Figure 23](#))



A **Select results folder** window appears.

2. Click to highlight the previously ran analysis folder you want to view in detail, then click **Select Folder**. **Note:** Past results can only be viewed if the original files (inside the folder you select) have not been modified.

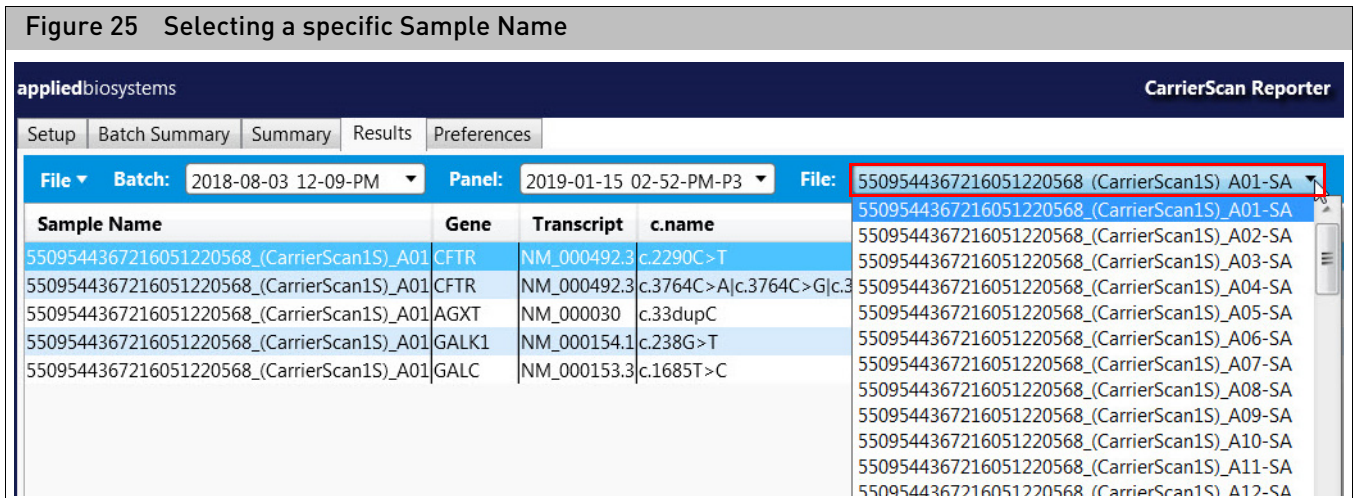
A detailed analysis of your selected folder appears in the Analysis Results Table, as shown in [Figure 24](#).

Figure 24 Results Table

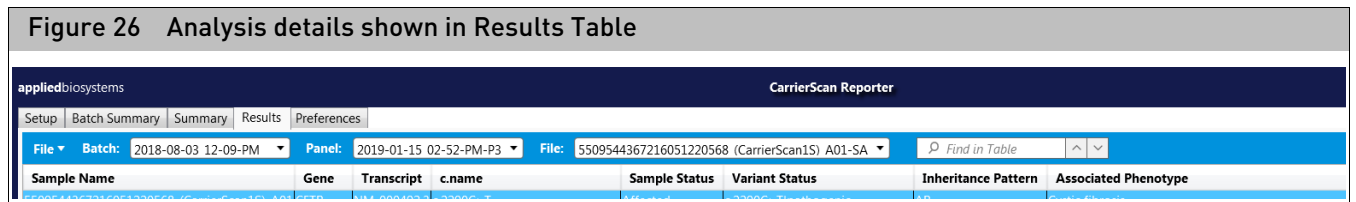
The screenshot shows the 'CarrierScan Reporter' software window with the 'Results' tab selected. The interface displays a table with the following data:

| Sample Name | Gene | Transcript | c.name | Sample Status | Variant Status | Inheritance Pattern | Associated Phenotype | Description | Severity | p.name |
|--|------|-------------|-------------------------------|---------------|----------------------|---------------------|--------------------------------|-------------|----------|------------------------|
| 5509544367216051220568_(CarrierScan1S)_A01 | CFTR | NM_000492.3 | c.2290C>T | Affected | c.2290C>Tpathogenic | AR | Cystic fibrosis | | B | p.Arg764Trp.Arg764Ter |
| 5509544367216051220568_(CarrierScan1S)_A01 | CFTR | NM_000492.3 | c.3764C>A/c.3764C>G/c.3764C>T | Carrier | c.3764C>A/pathogenic | AR | Cystic fibrosis | | B | p.Ser1255Leu.p.Ser1255 |
| 5509544367216051220568_(CarrierScan1S)_A01 | AGXT | NM_000030 | c.33dupC | Affected | c.33dupC/pathogenic | AR | Hyperoxaluria; primary; type 1 | | B | p.Lys12Glnfs |

- Click the **Sample Name** drop-down menu (Figure 25) to select a sample to view.



The Results table populates with your selected Sample Name’s details, as shown in Figure 26.



See the table below for column definitions.

| Column | Description |
|----------------------|--|
| Sample Name | Displays the filename of the sample. |
| Gene Name | Name of heritable genetic sequence that encodes proteins. |
| Transcript c.name | RefSeq transcript ID associated with the c.name. |
| Sample Status | Hereditary status of the sample for the associated phenotype based on information from the public domain. Note: An asterisk [*] indicates that the variant has varying calls and should be reviewed. |
| Variant Status | Standard classification of variant significance based on ACMG guidelines (pathogenic, likely pathogenic, etc.) Note: An asterisk [*] indicates variable or conflicting significance across reported sources. For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25741868 |
| Inheritance Pattern | Method of Inheritance (Example: AR (Autosomal Recessive), XLR (X-linked Recessive)) |
| Associated Phenotype | Displays the Phenotype that is associated with the variant. |
| Description | Detailed information about the Phenotype. |

| Column | Description |
|-----------------|---|
| Severity | Severity and impact rating of phenotypes. A = Profound, B = Severe, U = Unclassified, based on systematic classifications. For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25494330 |
| c.name | Displays standard variant nomenclature based on coding DNA reference sequences. |
| p.name | Displays the change in protein translation for the variant. |
| g.name | Displays the genomic coordinates for the variant. |
| Alternate Names | Displays the alternative names for the variant. |

Searching table content

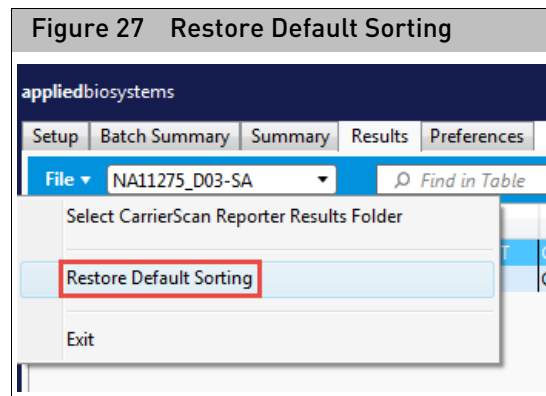
1. From the **Search** box in the Results window tab, type the name (or portion of the name) you want to locate.
2. Click the up or down arrow (or press the **Enter** key).
If a match is found, it is highlighted in the table.
3. Click the up or down arrow (or press the **Enter** key) to go to the next/previous matching entry.

Default row sorting

The Results table (Figure 26) displays Variant Statuses containing 'pathogenic' content first, then 'pseudo'-deficiency related variants.

Click once on a column header to sort its rows (in ascending or descending order) or drag-and-drop a row(s) to its new location within the table.

If you want to return to the default sorting order, click **File** → **Restore Default Sorting**.



Note: Custom sorting is not saved, as exiting and then reopening the application returns the Results table to its default sorted order.



Exporting

Exports

For each analysis run, four exports are auto-generated.

Each of these exports reside in the Output folder you assigned earlier [on page 8](#).

- "Full export"
- "Positive export" [on page 30](#)
- "Summary export" [on page 30](#)
- "Review export" [on page 31](#)

The exports for an Individual analysis provide information pertaining to the single sample in the analysis.

The exports for the Paired analysis provide information pertaining to both samples used in the paired analysis.

Full export

The Full Export details the data for all the markers in the panel whether or not a variant was detected. It also contains all probesets designed to analyze markers in question (especially for genotyping (GT) and Indel markers where multiple probesets are used to check for a particular SNP).

Note: The Full Export column definitions listed below were referenced from the White Paper *PMID: 25730230 DOI: 10.1097/AOG.000000000000666*. To access this publication, go to: <https://www.ncbi.nlm.nih.gov/pubmed/25730230>

IMPORTANT! The table below lists the columns included in the CarrierScan Reporter's commercial library file package. By default, all columns are displayed in all outputs and in the tables, but are not essential for the software to run. However, if you are creating a custom annotation panel, the columns noted with an asterisk [*] must be included in your annot.panel file.

| Column | Description |
|-------------------|--|
| Sample Name * | Displays the filename of the sample. |
| Gene Name * | Name of heritable genetic sequence that encodes proteins. |
| Transcript c.name | RefSeq transcript ID associated with the c.name. |
| Sample Status | Hereditary status of the sample for the associated phenotype based on information from the public domain. Note: An asterisk [*] indicates that the variant has varying calls and should be reviewed. |
| Variant Status * | Standard classification of variant significance based on ACMG guidelines (pathogenic, likely pathogenic, etc.) Note: An asterisk [*] indicates variable or conflicting significance across reported sources. For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25741868 |

| Column | Description |
|---|---|
| Inheritance Pattern | Method of Inheritance (Example: AR (Autosomal Recessive), XLR (X-linked Recessive)) |
| Associated Phenotype | Displays the Phenotype that is associated with the variant. |
| Description | Detailed information about the Phenotype. |
| Severity | Severity and impact rating of phenotypes based on these systematic classifications: A = Profound, B = Severe, U = Unclassified For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25494330 |
| c.name * | Displays standard variant nomenclature based on coding DNA reference sequences. |
| p.name | Displays the change in protein translation for the variant. |
| g.name | Displays the genomic coordinates for the variant. |
| Alternate Names | Displays the alternative names for the variant. |
| Carrier Frequency - Population 1 (Optional) | Proportion of the named Population 1 that has the recessive trait or phenotype. Note: This column is included if it is on the annotation panel used for analysis. |
| Detection Rate Population 1 (Optional) | Theoretical proportion of carriers in the named Population 1 that is identified by this assay. Note: This column is included if it is on the annotation panel used for analysis. |
| Residual Chance of Occurrence Population 1 (Optional) | Remaining chance the sample will be a carrier of the associated phenotype after a negative result considering Population 1's carrier frequency. For more information, see Appendix A, "Chance of occurrence calculation guidelines" . Note: This column is included if it is on the annotation panel used for analysis. |
| Marker Type | The type of marker (Indel, SNP, CN). |
| Affy SNP ID * | Unique Thermo Fisher Scientific generated identifier for the SNP. |
| Probeset ID * | Unique Thermo Fisher Scientific identifier for the probeset. |
| Recommended Probeset | A quality control metric determined by SNP Polisher algorithm that chooses the best probesets querying a SNP. |
| Probe Count | The number of probes used for the call. |
| Confidence | The number of probesets that agree with determined sample status for that variant. |
| Analysis Call | Call code for the variant (AA, AB, BB, etc). |
| Translated Genotype Call | The nucleotide conversion of the genotyping call code. |
| Ref Code | The Call Code (A,B,C) of the Reference Allele. |
| Alt Code | The Call Code (A,B,C,D,.....) of the Alternate Allele. |
| Ref Allele * | The call for the reference allele associated with a normal phenotype. |
| Alt Allele * | The call for the 1st alternate allele associated with a non-normal phenotype. |
| Variant Status Alt Allele | Severity status for the variant mapped to Alt Allele. |
| CN Region | The defined region used for copy number analysis. |
| CN State | The copy number state of the defined region. |
| Chromosome * | The chromosomal location of the variant. |
| Physical Position * | Starting genomic coordinates of the variant bases. |

| Column | Description |
|---------------|--|
| Position End | Ending genomic coordinates of the variant bases. |
| Plate Barcode | Barcode of the array plate. |
| Well Position | Position location of the array on the plate. |

Positive export

Positive Export details the data for ALL markers in the panel that have a non-normal genetic state (positive result).

Note: Positive Export is a subset of the Full Export, therefore refer to the table of definitions above, as the columns featured in the Full Export and Positive Export are identical.

Summary export

The Summary Export provides summary results for ALL phenotypes in the panel that have at least one non-normal genetic event, including Affected, Carrier, and CNgain sample statuses associated with pathogenic and pseudoallele/pseudodeficiency related variant statuses.

The Summary Export contains the information provided in the Positive Export collated by variant. This export is also read into the software and shown in the Analysis Results Table.

The Positive Export and Summary Export column definitions are listed in the table below.

| Column | Description |
|----------------------|--|
| Sample Name | Displays the filename of the sample. |
| Gene Name | Name of heritable genetic sequence that encodes proteins. |
| Transcript c.name | RefSeq transcript ID associated with the c.name. |
| Variant Status | Standard classification of variant significance based on ACMG guidelines (pathogenic, likely pathogenic, etc.) Note: An asterisk [*] indicates variable or conflicting significance across reported sources. For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25741868 |
| Sample Status | Hereditary status of the sample for the associated phenotype based on information from the public domain. Note: An asterisk [*] indicates that the variant has varying calls and should be reviewed. |
| Inheritance Pattern | Method of Inheritance (Example: AR (Autosomal Recessive), XLR (X-linked Recessive)) |
| Associated Phenotype | Displays the Phenotype that is associated with the variant. |
| Description | Detailed information about the Phenotype. |
| Severity | Severity and impact rating of phenotypes. A = Profound, B = Severe, U = Unclassified, based on systematic classifications. For more information, go to: https://www.ncbi.nlm.nih.gov/pubmed/25494330 |
| c.name | Displays standard variant nomenclature based on coding DNA reference sequences. |
| p.name | Displays the change in protein translation for the variant. |
| g.name | Displays the genomic coordinates for the variant. |

| Column | Description |
|-----------------|---|
| Alternate Names | Displays the alternative names for the variant. |
| Plate Barcode | Barcode of the array plate. |
| Well Position | Position location of the array on the plate. |

Review export

The Review Export provides details for all variants that did not have a result on the panel. If a variant has a Sample Status result of Normal, Carrier, Affected, CN gain assessment, then it is considered tested. Otherwise, the variant is recorded in this report to be reviewed.

The Review Export column definitions are listed below.

| Column | Description |
|---------------|---|
| Sample Name | Displays the filename of the sample. |
| Gene Name | Name of heritable genetic sequence that encodes proteins. |
| c.name | Displays standard variant nomenclature based on coding DNA reference sequences. |
| p.name | Displays the change in protein translation for the variant. |
| Probeset ID | Unique Thermo Fisher Scientific identifier for the probeset. |
| CN Region | The defined region used for copy number analysis. |
| Failure Mode | Reason the Probeset or CN Region failed to give a sample status. |
| Plate Barcode | Barcode of the array plate. |
| Well Position | Position location of the array on the plate. |



Chance of occurrence calculation guidelines

Single and paired sample analysis reports contain ethnic panels pre-populated with known frequencies and detection rates.

Note: The Chance of Occurrence Guideline calculations (shown below) were referenced from the European Journal of Human Genetics White Paper *PMID: 18685558* *PMCID: PMC2985951* *DOI: 10.1038/ejhg.2008.136*. To access this publication, go to: <https://www.ncbi.nlm.nih.gov/pubmed/18685558>

- Calculations for paired residual chance of occurrence (RCO) are:
 $(RCO1) \times (RCO2) \times 1/4 = \text{paired RCO}$
 - **RCO1** = Residual chance of occurrence for sample 1 for the assigned ethnicity and associated phenotype of normal sample status.
 - **RCO2** = Residual chance of occurrence for sample 2 for the assigned ethnicity and associated phenotype of normal sample status.
 - When the sample status is deemed **Carrier** or **Affected**, the values used for RCO are **1** and **2** respectively.

Special handling is performed for Alpha-Thalassemia sample status determination that affects the RCO values used in calculations. Copy number states of HBA1 and HBA2 are considered together (according to guidelines referenced from NCBI GeneReviews). To access this publication, go to: <https://www.ncbi.nlm.nih.gov/books/NBK1435>

For example, in all reports (exclusive of the BatchSummaryResults) the status for both HBA1 and HBA2 copy number regions will always have the same sample status regardless if the genetic aberration occurs on only 1 region. In the BatchSummaryResults file, the HBA1 and HBA2 copy number regions are assessed independently to allow for review of each variant.

In the scenario of a cis/trans single deletion of HBA1 and HBA2, the cis orientation is assumed as input into the paired RCO calculations.

For support visit [thermofisher.com/support](https://www.thermofisher.com/support) or email techsupport@lifetech.com
thermofisher.com

4 June 2019

ThermoFisher
SCIENTIFIC