

Genotyping Analysis Module

USER GUIDE

for use with QuantStudio™ Design and Analysis Software v2

Publication Number MAN0018749

Revision D.0



Revision history: MAN0018749 D.0 (English)

Revision	Date	Description
D.0	15 December 2023	<p>The document was updated for QuantStudio™ Design and Analysis Software v2.8 and the Genotyping Analysis Module v1.6.</p> <ul style="list-style-type: none">• Permissions were added for the following tasks, when the security settings are enabled:<ul style="list-style-type: none">– Selecting the location when a file is saved (“Select a system template or existing plate file to set up a new plate file” on page 7).– Selecting a passive reference (“Select a passive reference” on page 11).• Instructions were added to export the results (“Export the results” on page 16).
C.0	5 July 2023	<p>The document was updated for QuantStudio™ Design and Analysis Software v2.7 and the Genotyping Analysis Module v1.5.</p> <ul style="list-style-type: none">• The list of compatible data files was removed. See the documentation for the main software for this information.• Information was added about the OpenArray™ Plate format.• The instructions to select the genotyping analysis module were updated to indicate that OpenArray™ Plate genotyping data files automatically have the genotyping analysis module applied when they are opened.
B.0	15 April 2020	Remove send to the instrument run queue.
A.0	14 October 2019	New document.

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.

NOTICE TO PURCHASER: DISCLAIMER OF LICENSE: Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Thermo Fisher Scientific, either expressly, or by estoppel.

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a trademark of Roche Molecular Systems, Inc., used under permission and license.

©2019-2023 Thermo Fisher Scientific Inc. All rights reserved.

Contents

- **CHAPTER 1** About the Genotyping Analysis Module 5
- **CHAPTER 2** Workflow: Genotyping analysis 6
- **CHAPTER 3** Set up a plate file for genotyping analysis 7
 - Select a system template or existing plate file to set up a new plate file 7
 - Confirm or edit the run method for genotyping analysis 8
 - Confirm or edit the plate setup for genotyping analysis 8
 - Add samples and assign to wells 9
 - Add SNP assays and assign to wells 9
 - Edit reagent information 10
 - Select a passive reference 11
 - Review and save the plate file 11
- **CHAPTER 4** Perform genotyping analysis 12
 - Review results in the **Amplification Plot** 12
 - Select the Genotyping Analysis Module 12
 - Review results in the **Allelic Discrimination Plot** 12
 - Omit outliers from genotyping analysis 14
 - (Optional)* Review dye signal profile in the **Multicomponent Plot** 15
 - (Optional)* Review signal profile in the **Raw Data Plot** 15
 - Edit **Genotyping Analysis Setting** 15
 - Export the results 16
- **CHAPTER 5** About genotyping analysis 17
 - Overview of genotyping analysis 17
 - Sample types for genotyping analysis 17
 - Allelic Discrimination Plot overview 18

- **APPENDIX A Documentation and support** 19
 - Related documentation 19
 - Customer and technical support 19
 - Limited product warranty 19
- Index** 20



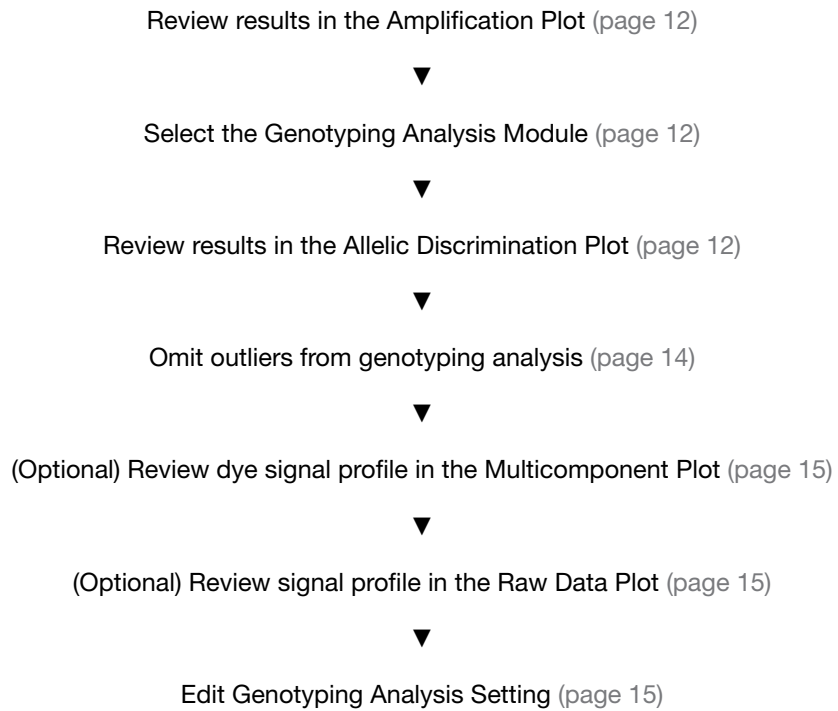
About the Genotyping Analysis Module

The Genotyping Analysis Module for QuantStudio™ Design and Analysis Software v2 is used to detect single nucleotide polymorphism (SNP) variants of a target nucleic acid sequence.

For more information about genotyping analysis, see Chapter 5, “About genotyping analysis”.

2

Workflow: Genotyping analysis



IMPORTANT! If you omit wells or configure the analysis settings, click **Analyze** to reanalyze the data.



Set up a plate file for genotyping analysis

For detailed instructions about setting up a plate file, see [? Help ▶ Help Contents](#).


Select a system template or existing plate file to set up a new plate file

A plate file contains the information that is necessary to perform an instrument run, including instrument setup, run method, plate setup, and analysis setting.

A system template is a non-editable plate file that is included with the software.

A new plate file must be created from a system template or a previously created plate file.

For detailed information about system templates and plate files, see [? Help ▶ Help Contents](#).

1. In the home screen, click  **Set Up Plate**.
The **Plate Gallery** opens to the **System Templates** tab.

2. **IMPORTANT!** Select a system template or a plate file that corresponds to your instrument, block, and run mode. These properties are not editable after the plate file has been created.

In the left pane, select the appropriate options to filter the system template and plate file lists.

- **Instrument**
- **Block**
- **Run Mode**
- **Analysis**

Note: Thermal protocol, plate setup, and post-run analysis options are independent of analysis module selection. Analysis module selection can be changed at any point during plate file set or post-run analysis (see “Select the Genotyping Analysis Module” on page 12).

3. Navigate to, then select a system template or plate file.

Tab	Description
System Templates	Contains system templates, non-editable plate files that are included with the software. Select a system template to automatically generate a new plate file that can be edited, then saved.
My Plate Files	Contains plate files that were previously saved to My Plate Files . plate files that are included with the software. Select an existing plate file to edit, then save, or to save as a new plate file.
Recents	Contains plate files that were recently opened. Recently opened plate files from System Templates and My Plate Files do not populate this tab. Select an existing plate file to edit, then save, or to save as a new plate file.

Note: To view all options for opening the plate file, hover over the plate file, then click ... **(Actions)**.

Selecting the location where the file is saved is a controlled function. If the **Browse** button in the **Save As** dialog box is inactive, the security has been enabled in the software. The user role does not have the permission to select the location where the file is saved. The user role must have the permission of **Edit Save As Destination**.

The plate file opens in the **Run Method** tab.

Confirm or edit the run method for genotyping analysis

For most analysis, the default run method is appropriate. The following options are compatible for genotyping analysis.

- PCR
- 1-step RT-PCR
- 2-step RT-PCR
- In a plate file, in the **Run Method** tab, adjust the run method elements as needed.
For detailed instructions about editing the run method, see [Help](#) ▶ **Help Contents**.
- Click ... **(Actions)** ▶ **Filter Settings** to confirm or edit filter settings.

Confirm or edit the plate setup for genotyping analysis

For detailed instructions about plate setup, or to download example plate setup files, see [Help](#) ▶ **Help Contents**.

Add samples and assign to wells

For detailed instructions about plate setup, see [?](#) **Help** ▶ **Help Contents**.

1. In the **Plate Setup** tab, add samples and assign to wells using the following options.

- Import a plate setup file
- Manually add samples to the **Samples** table
- Manually add samples to wells in the plate layout

For the OpenArray™ Plate format, a sample layout can be imported. Samples cannot be added manually. The other formats allow a plate setup file to be imported, including targets and assays.

2. Confirm or edit sample information in the **Samples** table.

Column	Description
Name	Sample name
Color	Sample color
Type ^[1]	Genotyping analysis uses the following sample types. <ul style="list-style-type: none"> • Unknown • Positive 1/1 • Positive 1/2 • Positive 2/2 • Negative Control

^[1] For more information, see “Sample types for genotyping analysis” on page 17.

3. Confirm or edit sample well assignments in the plate layout.

Add SNP assays and assign to wells

For detailed instructions about plate setup, see [?](#) **Help** ▶ **Help Contents**.

SNP assays cannot be imported or assigned for the OpenArray™ Plate format.

1. In the upper-left corner of the **Plate Setup** tab, click **SNP**.

2. Add SNP assays and assign to wells using the following options:

- Import an Assay Information File (AIF)
- Import a plate setup file
- Manually add SNP assays to the **SNP Assays** table
- Manually add SNP assays to wells in the plate layout
- Import TaqMan™ assay plate and card files

3. Confirm or edit target information in the **SNP Assays** table.

Column	Description
Name	SNP assay name
Color	SNP assay color
Allele 1 Reporter	Fluorescent reporter dye for allele 1 of the SNP assay
Allele 1 Quencher	Quencher for allele 1 of the SNP assay
Allele 2 Reporter	Fluorescent reporter dye for allele 2 of the SNP assay
Allele 2 Quencher	Quencher for allele 2 of the SNP assay
Task ^[1]	<p>The software automatically assigns a task to the target in a well based on the sample type in that well. The following tasks are used for standard curve analysis.</p> <ul style="list-style-type: none"> • Unknown • Positive 1/1 • Positive 1/2 • Positive 2/2 • Negative Control

^[1] For more information, see "Sample types for genotyping analysis" on page 17.

4. Confirm or edit SNP assay well assignments in the plate layout.

Edit reagent information

Reagents can be edited for the OpenArray™ Plate format.

1. In the **Plate Setup** tab, in the **Targets/SNP Assays** table pane, click **Reagents**.
2. In the **Reagents** table, click **+ (Add)**.

Note:

- Click **⋮ (Actions) ▶ Export Reagents** to export reagents.
 - Click **⋮ (Actions) ▶ Import Reagents** to import reagents.
 - Click **⋮ (Actions) ▶ Scan Reagents** to scan reagents.
-

3. Enter the following information for each reagent.

- | | |
|-----------|-------------------|
| • Name | • Part Number |
| • Type | • Lot Number |
| • Barcode | • Expiration Date |
-

Note: If the master mix that you enter is not compatible with the current run method, you have the option to apply the recommended run method for your master mix, instrument, block, and run mode.

4. (Optional) Click **✕ (Remove)** in the row of a reagent to delete it from the table.

Select a passive reference

Editing the passive reference is a controlled function. If security has been enabled in the software, the user role must have the permission of **Edit Passive Reference**.

1. In the upper-left corner of the **Plate Setup** tab, select a passive reference from the dropdown list.
2. *(Optional)* Save the plate file or data file.

Review and save the plate file

1. In the **Run Summary** tab, review the run method selections, then edit if needed.
2. Review the plate setup, then edit if needed.
3. *(Optional)* Click the barcode field, then scan the plate barcode.
4. *(Optional)* Select **Add to My Plates**.
This option allows you to create new plate files using the current plate file as a template.
5. Select an instrument from the list.
If the instrument does not appear on the list, click **System ▶ Instruments** to add a new instrument.
6. Save the plate file.

Start the run on an instrument. For specifics on starting an instrument run, see the instrument documentation.

4

Perform genotyping analysis

Review results in the Amplification Plot

For detailed instructions about reviewing results in the **Amplification Plot**, see [?](#) **Help** ▶ **Help Contents**.

If no data are displayed in the **Quality Check** tab, or if reanalysis is required, click **Analyze**.

1. In the **Quality Check** tab, in the plot pane, select **Amplification Plot** from the dropdown list.
2. Review the amplification status for each well.
3. Review or edit threshold settings.
4. Review or edit baseline settings.

Select the Genotyping Analysis Module

OpenArray™ Plate genotyping data files automatically have the genotyping analysis module applied when they are opened.

1. In an open data file, click **Actions** ▶ **Analysis Modules**.
2. In the **Analysis Modules** window, select **Genotyping**, then click **Ok**.
The Genotyping Analysis Module opens.

Click **Analyze**, then review the results in the **Genotyping** tab.

Review results in the Allelic Discrimination Plot

For more information about the allelic discrimination plot, see “Allelic Discrimination Plot overview” on page 18.

If no data are displayed in the **Genotyping** tab, or if reanalysis is required, click **Analyze**.

1. In the **Genotyping** tab, in the upper-left corner of the **Allelic Discrimination Plot** pane, select an assay from the **Assays** list (see Figure 1 on page 14).
The Allelic Discrimination Plot is displayed for the selected assay.

2. Confirm that the data points for each genotype control cluster along the expected axis of the plot.

Table 1 Cluster assignments in an allelic discrimination plot

Content of samples	Location in AD plot
Allele 1 homozygote ^[1]	Lower right corner
Allele 2 homozygote ^[2]	Upper left corner
Alleles 1/2 heterozygote ^[3]	Approximately midway between Allele 1 and Allele 2 clusters
No-template control	Bottom left corner
Undetermined	Anywhere outside the regions described above
No amplification	With no-template control (NTC) cluster in the bottom left corner

^[1] Labeled with allele 1 fluorescent dye.

^[2] Labeled with allele 2 fluorescent dye.

^[3] Labeled with allele 1 and allele 2 fluorescent dyes.

3. Confirm that the cluster at the bottom-left corner of the plot contains only negative control wells. Samples can unexpectedly cluster with the negative controls for one of the following reasons.
 - Samples contain no DNA
 - Samples contain PCR inhibitors
 - Samples are homozygous for a sequence deletion
4. Visually evaluate clusters for the three possible genotypes (see Figure 1 on page 14).

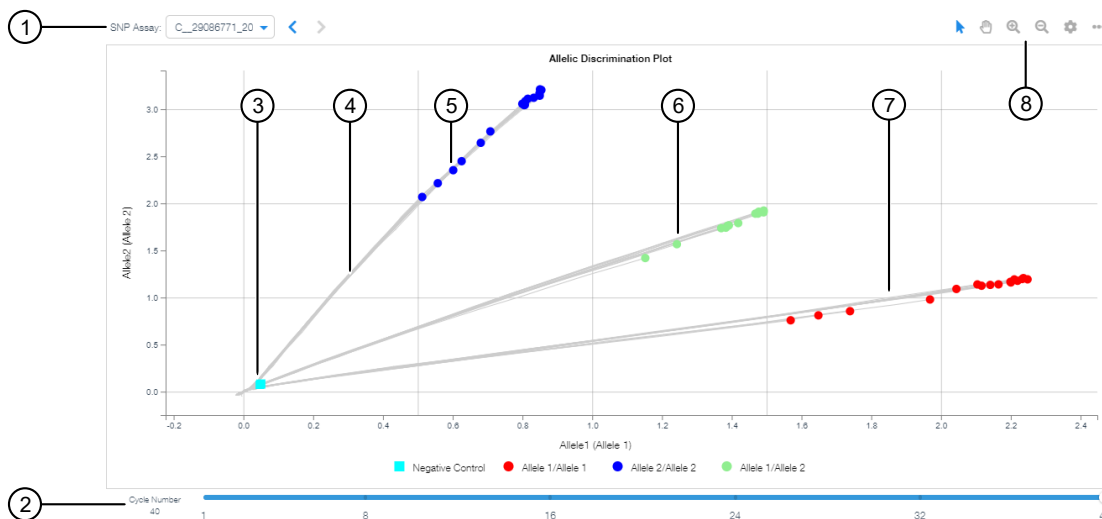


Figure 1 Allelic Discrimination Plot

- ① Assay selection
- ② **Cycle Number** slider—For analysis that includes real-time data, the slider allows you to display the allelic discrimination data for all samples at each cycle during the PCR amplification.
- ③ NTC cluster
- ④ Real-time PCR amplification trails—For analysis that includes real-time data
- ⑤ Allele 2 homozygote cluster
- ⑥ Allele 1/2 heterozygote cluster
- ⑦ Allele 1 homozygote cluster
- ⑧ Plot toolbar

To confirm results, retest outliers and samples with no amplification (cluster with negative controls).

Omit outliers from genotyping analysis

Outlier wells have C_q values that differ significantly from the average for the associated replicate wells. To ensure C_q precision, consider omitting the outliers from analysis.

1. In the **Genotyping** tab, select an option to omit wells from analysis.

Option	Description
Omit wells in the Plate Layout	Select outlier wells, then click ⋮ (Actions) ▶ Omit Wells .
Omit wells in the Well Table	Select Omit in the row of the outlier well.

2. Click **Analyze** to reanalyze the data with any outliers removed.

(Optional) Review dye signal profile in the Multicomponent Plot

For more information about the **Multicomponent Plot**, see [?](#) **Help** ▶ **Help Contents**.

If no data are displayed in the **Quality Check** tab, or if reanalysis is required, click **Analyze**.

1. In the **Quality Check** tab, in the plot pane, select **Multicomponent Plot** from the dropdown list.
2. Review the signal profiles for the passive reference dye, reporter dye, and negative control wells.
3. Review the plot to ensure that there are no irregularities in the dye signals.

(Optional) Review signal profile in the Raw Data Plot

For detailed instructions about reviewing results in the **Raw Data Plot**, see [?](#) **Help** ▶ **Help Contents**.

If no data are displayed in the **Quality Check** tab, or if reanalysis is required, click **Analyze**.

1. In the **Quality Check** tab, in the plot pane, select **Raw Data Plot** from the dropdown list.
2. Click-drag the **Cycle Number** slider through all of the cycles, then confirm that each filter displays the characteristic signal increase.

Edit Genotyping Analysis Setting

Open the Genotyping Analysis Module.

1. Click **Actions** ▶ **Genotyping Analysis Setting**.
2. Select an option from the **Analysis Data** list to determine the data that used for genotyping analysis.

Analysis Data	Description
Post Read	Only post-PCR read data is used to determine calls.
Pre Read & Post Read ^[1]	The pre-PCR read is subtracted from the post-PCR read to determine calls.
Real Time Data ^[2]	The normalized reporter data (Rn) from the selected cycle of the cycling stage is used to determine calls. Analysis is not restricted to the last cycle. Adjust the analysis cycle using the Cycle Slider in the Allelic Discrimination Plot pane (see Figure 1 on page 14).

^[1] The run method must include a pre-read stage.

^[2] Data collection must be on during the PCR stage.

3. (Optional) Edit the following settings:

- **Min Confidence %**: Change the confidence threshold to increase or decrease stringency for making a genotype call.
- (Real Time Data Analysis Only) **Baseline Start** and **Baseline End**: Enter the first and last cycle that the software will use to calculate the baseline.
- (Real Time Data Analysis Only) **Cycle Number**: Change the cycle number used to generate the normalized reporter data (Rn).

Option	Description
Edit default analysis settings	<ol style="list-style-type: none"> 1. In the table, in the Default Setting row, click in a cell to edit the setting. 2. To apply the default analysis settings to an assay, in the assay row, select Use Default.
Edit analysis settings for an individual assay	In the table, in the assay row, deselect the Use Default checkbox, then click in a cell to edit the setting.

4. Click **Apply**.

The data is reanalyzed using the updated analysis settings.

Export the results

1. In the table pane, click **⋮ (Actions) ▶ Export**.
2. (Optional) In the **Export CSV** dialog box, edit the file name in the **File Name** field.
The **File Name** field is populated with a default file name.
3. Click **Browse** to select a location to save the file.
Selecting the location where the file is saved is a controlled function. If the **Browse** button in the **Export CSV** dialog box is inactive, the security has been enabled in the software. The user role does not have the permission to select the location where the file is saved. The user role must have the permission of **Edit Export Destination**.
4. Click **Export**.



About genotyping analysis

Overview of genotyping analysis

Genotyping analysis is used to detect single nucleotide polymorphism (SNP) variants of a target nucleic acid sequence.

Genotyping analysis uses preformulated TaqMan™ SNP Genotyping Assays that include the following components:

- Two sequence-specific primers for amplification of sequences containing the SNP of interest
- Two allele-specific TaqMan™ probes for Allele 1 and Allele 2

For genotyping analysis, the software performs the following tasks:

1. The software normalizes the fluorescence of the reporter dyes to the fluorescence of the passive reference dye in each well.
2. The software plots the normalized reporter dye signal of each sample well on an Allelic Discrimination Plot, which contrasts the reporter dye intensities of the allele-specific probes.
3. The software algorithmically clusters the sample data, and assigns a genotype call to the samples of each cluster according to its position on the plot.

Sample types for genotyping analysis

Genotyping analysis includes the following sample types.

Sample type (Type column in Samples table)	Sample description	Automatic target task assignment ^[1] (Task column in Targets table)
Unknown	Test sample	Unknown
Positive 1/1	Control sample that is heterozygous allele 1/allele 2	Positive 1/1
Positive 1/2	Control sample that is heterozygous allele 1/allele 2	Positive 1/2
Positive 2/2	Control sample that is homozygous for allele 2	Positive 2/2
Negative Control	Water or buffer No amplification of the target should occur in NTC wells.	Negative control

^[1] The software automatically assigns a task to the target in a well based on the sample type in that well.

Allelic Discrimination Plot overview

The Allelic Discrimination Plot contrasts the R_n or the ΔR_n of the reporter dyes for the allele-specific probes of the SNP assay. It is an intermediary step in the software algorithm for genotyping calls.

A typical Allelic Discrimination Plot shows homozygote clusters, a heterozygote cluster, and the no-template controls. The points in each cluster are grouped closely together and each cluster is located well away from the other clusters. For expected cluster locations in the plot, see Figure 2 on page 18.

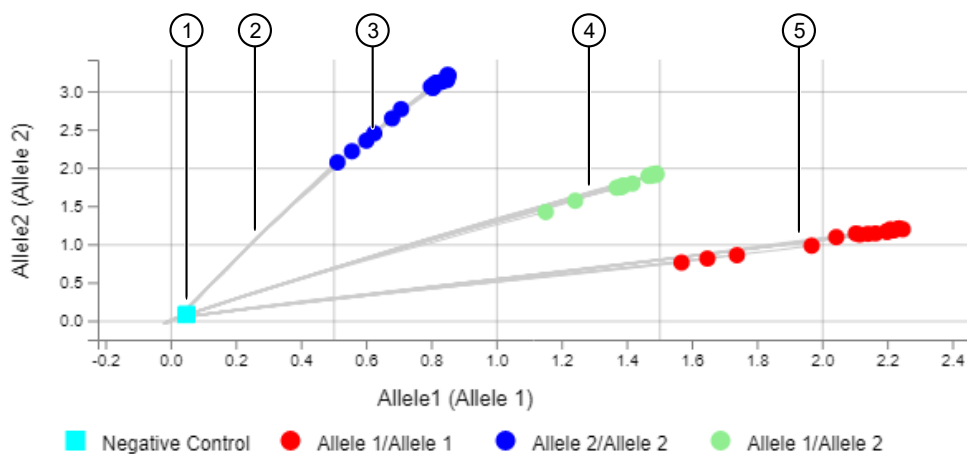


Figure 2 Example allelic discrimination plot

- ① No-template control cluster
- ② Real-time PCR amplification trails (for analysis that includes real-time data)
- ③ Allele 2 homozygote cluster
- ④ Allele 1/2 heterozygote cluster
- ⑤ Allele 1 homozygote cluster

Note:

- Undetermined samples may appear anywhere on the plot.
 - No-amplification samples cluster with the no-template control samples.
-

Note: The clustering algorithm does not call genotypes if all the samples are one genotype (form one cluster).



Documentation and support

Related documentation

Document	Publication number
<i>QuantStudio™ Design and Analysis Software v2 User Guide</i>	MAN0018200
<i>QuantStudio™ Design and Analysis v2 User Guide (Thermo Fisher™ Connect Platform)</i>	MAN0018202

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation, including:
 - 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 found at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at <http://www.thermofisher.com/support>.

Index

A

Allelic Discrimination Plot
 overview 18
 review results 12
Amplification Plot, review 12
analysis module, select 12

G

genotyping analysis
 overview 17
 perform 12
 sample types 17
 samples 9
 SNP assays 9
 workflow 6
genotyping analysis setting, edit 15
genotyping experiments, Allelic Discrimination Plot
 18

L

limited product warranty 19

M

MultiComponent Plot, review 15

O

omit wells, genotyping analysis 14

P

passive reference, select 11

plate file
 analysis type 7
 create 7
 save 11
 set up 7, 8
plate setup
 samples 9
 SNP assays 9

R

Raw Data Plot, review 15
reagents, information 10
related documentation 19
run method 8

S

sample types, genotyping analysis 17
samples 9
SNP assays 9
support, customer and technical 19

T

terms and conditions 19

W

warranty 19
workflow 6

