

# Standard Curve Analysis Module

## USER GUIDE

for use with:

QuantStudio™ Design and Analysis Software v2

Publication Number MAN0018746

Revision B.0



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

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.

**Revision history:** Pub. No. MAN0018746

Revision	Date	Description
B.0	15 April 2020	Changes for version 1.4: Remove send to the instrument run queue; add Replicate Group Table.
A.0	26 August 2019	New document.

**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 registered trademark of Roche Molecular Systems, Inc., used under permission and license.

©2020 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

■	<b>CHAPTER 1</b>	About the Standard Curve Analysis Module .....	5
		Compatible data files .....	5
■	<b>CHAPTER 2</b>	Workflow: Standard curve analysis .....	7
■	<b>CHAPTER 3</b>	Set up a plate file for standard curve analysis .....	8
		Select a system template or existing plate file to set up a new plate file .....	8
		Confirm or edit the run method for standard curve analysis .....	9
		Confirm or edit the plate setup for standard curve analysis .....	9
		Add samples and assign to wells .....	10
		Add targets and assign to wells .....	10
		Set up the standard curve .....	11
		Edit reagent information .....	12
		Select a passive reference .....	12
		Review and save the plate file .....	12
■	<b>CHAPTER 4</b>	Perform standard curve analysis .....	14
		Review results in the <b>Amplification Plot</b> .....	14
		Select the Standard Curve Analysis Module .....	14
		Review results in the <b>Standard Curve Plot</b> .....	14
		Identify and omit outliers from standard curve analysis .....	15
		(Optional) Review dye signal profile in the <b>MultiComponent Plot</b> .....	15
		(Optional) Review signal profile in the <b>Raw Data Plot</b> .....	16
		(Optional) Edit standard curve analysis settings .....	16
■	<b>CHAPTER 5</b>	About standard curve analysis .....	17
		Overview of standard curve analysis .....	17
		Sample types for standard curve analysis .....	17
		<b>Standard Curve Plot</b> overview .....	18

Documentation and support .....	20
Related documentation .....	20
Customer and technical support .....	20
Limited product warranty .....	20
Index .....	21



# About the Standard Curve Analysis Module

The Standard Curve Analysis Module for QuantStudio™ Design and Analysis Software v2 is used to determine absolute target quantity in test samples.

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

## Compatible data files

The software is compatible with data files for the following real-time PCR instruments:

- QuantStudio™ 7 Pro Real-Time PCR System
- QuantStudio™ 6 Pro Real-Time PCR System
- QuantStudio™ 5 Real-Time PCR System (if the plate file for the instrument run is created using QuantStudio™ Design and Analysis Software v2)
- QuantStudio™ 3 Real-Time PCR System (if the plate file for the instrument run is created using QuantStudio™ Design and Analysis Software v2)
- QuantStudio™ 1 Real-Time PCR System (if the plate file for the instrument run is created using QuantStudio™ Design and Analysis Software v2)

Data files for the following instruments are a legacy file format that can be opened in the software, but can only be saved as the updated file format:

- QuantStudio™ 5 Real-Time PCR System (if the plate file for the instrument run was not created using QuantStudio™ Design and Analysis Software v2)
- QuantStudio™ 3 Real-Time PCR System (if the plate file for the instrument run was not created using QuantStudio™ Design and Analysis Software v2)
- QuantStudio™ 1 Real-Time PCR System (if the plate file for the instrument run was not created using QuantStudio™ Design and Analysis Software v2)
- QuantStudio™ 6 Flex Real-Time PCR System
- QuantStudio™ 7 Flex Real-Time PCR System
- QuantStudio™ 12K Flex Real-Time PCR System (except OpenArray™ format)
- StepOnePlus™ Real-Time PCR System
- ViiA™ 7 Real-Time PCR System
- 7500/7500 Fast Real-Time PCR System
- 7900HT Real-Time PCR System



**Note:** To convert a legacy data file into the updated file format, open the data file, then click **Actions** ▶ **Save As**.



# Workflow: Standard curve analysis

For detailed instructions about setting up a plate file, or reviewing data in the **Quality Check** tab, see [?](#) **Help** ▶ **Help Contents**.

## Set up a plate file for standard curve analysis (page 8)

Select a system template or existing plate file to set up a new plate file (page 8)



Confirm or edit the run method for standard curve analysis (page 9)



Confirm or edit the plate setup for standard curve analysis (page 9)



Review and save the plate file (page 12)



## Perform standard curve analysis (page 14)

Review results in the Amplification Plot (page 14)



Select the Standard Curve Analysis Module (page 14)



Review results in the Standard Curve Plot (page 14)



Identify and omit outliers from standard curve analysis (page 15)



(Optional) Review dye signal profile in the MultiComponent Plot (page 15)



(Optional) Review signal profile in the Raw Data Plot (page 16)



(Optional) Edit standard curve analysis settings (page 16)



# Set up a plate file for standard curve 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 Standard Curve Analysis Module” on page 14).



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

Tab	Description
<b>System Templates</b>	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.
<b>My Plate Files</b>	Contains plate files that were previously saved to <b>My Plate Files</b> . 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.
<b>Recents</b>	Contains plate files that were recently opened. Recently opened plate files from <b>System Templates</b> and <b>My Plate Files</b> 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, mouse over the plate file, then click ... **(Actions)**.

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

## Confirm or edit the run method for standard curve analysis

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

**Note:** Multiple targets can be assayed using standard curve analysis, but each target requires its own standard curve. To set up the standard curve using the **Standard Curve Wizard** see “Set up the standard curve” on page 11.

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

Column	Description
Name	Sample name
Color	Sample color
Type <sup>[1]</sup>	Standard curve analysis uses the following sample types. <ul style="list-style-type: none"> <li>• Standard<sup>[2]</sup></li> </ul> <p><b>Note:</b> You must enter the quantity for each standard sample in the <b>Quantity</b> column.</p> <ul style="list-style-type: none"> <li>• Unknown</li> <li>• Negative Control</li> </ul>
Quantity (standard samples only)	Enter the quantity for the standard sample. <p><b>Note:</b> The quantity entered for a standard sample in the <b>Samples</b> table is used to populate the <b>Quantity</b> column for standard tasks in the <b>Targets</b> table (see “Add targets and assign to wells” on page 10).</p>

<sup>[1]</sup> For more information, see “Sample types for standard curve analysis” on page 17.

<sup>[2]</sup> Each target requires its own standard curve.

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

## Add targets and assign to wells

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

**Note:** Multiple targets can be assayed using standard curve analysis, but each target requires its own standard curve. To set up the standard curve using the **Standard Curve Wizard** see “Set up the standard curve” on page 11.

1. In the **Plate Setup** tab, add targets and assign to wells using the following options.
  - Import an AIF file
  - Import a plate setup file
  - Manually add targets to the **Targets** table
  - Manually add targets to wells in the plate layout

- Import TaqMan™ assay plate and card files

2. Confirm or edit target information in the **Target** table.

Column	Description
Name	Target name
Color	Target color
Task <sup>[1]</sup>	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. <ul style="list-style-type: none"> <li>• Standard</li> <li>• Unknown</li> <li>• Negative Control</li> </ul>
Quantity (standard tasks only)	The quantity entered for a standard sample in the <b>Samples</b> table is used to populate the <b>Quantity</b> column for standard tasks in the <b>Targets</b> table.

<sup>[1]</sup> For more information, see “Sample types for standard curve analysis” on page 17.

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

## Set up the standard curve

**Note:**

- Multiple targets can be assayed using standard curve analysis, but each target requires its own standard curve.
- To import a standard curve from a different data file, see “(Optional) Edit standard curve analysis settings” on page 16.
- You can also set up the standard curve during sample setup (see “Add samples and assign to wells” on page 10).

1. In the **Plate Setup** tab, in the plate setup pane, click ... **(Actions)** ▶ **Standard Curve Setup**.

The **Standard Curve Wizard** opens.

2. In the **Standard Curve Wizard** pane, enter the sample name prefix.

3. Select the target for the standard curve.

Option	Instructions
Target previously defined	Select the target from the dropdown list.
Target not previously defined	<ol style="list-style-type: none"> <li>1. Type the target name, the press <b>Enter</b>.</li> <li>2. Select a reporter from the dropdown list.</li> <li>3. Select a quencher from the dropdown list.</li> </ol>

4. Adjust the parameters for the dilution series if needed.
  - **Number of points**—5 recommended
  - **Number of replicates**—3 recommended
  - **Starting Quantity**—The highest or lowest standard quantity, without units.  
**Note:** The quantity must be  $\geq 1$ .
  - **Serial Factor**  
**Note:** The serial factor calculates quantities for all standard curve points.
    - Starting quantity is the highest value—Select 1:10 to 1:2.
    - Starting quantity is the lowest value—Select 2x to 10x.
5. Select an option to select the wells for the standard
  - Select **Automatically**.
  - Select **Manually**, then select wells using the displayed plate layout.
6. Select to arrange the standards in **Rows** or **Columns**.
7. Click **Apply Standard Curve**, then click **Close** to return to the **Plate Setup** tab.

## Edit reagent information

1. In the **Plate Setup** tab, in the **Targets/SNP Assays** table pane, click **Reagents**.
2. In the **Reagents** table, click **+** (**Add**).
3. Enter the reagent type, name, part number, lot number, and 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

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**.

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 standard curve 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 Standard Curve Analysis Module

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

Click **Analyze**, then review the results in the **Standard Curve** tab.

### Review results in the Standard Curve Plot

1. In the **Standard Curve** tab, in the plot pane, select a target from the **Targets** dropdown list.
2. In the plot pane, click , then select an option from the **Color By** dropdown list: **Target**, **Sample**, or **Task**.  
The plot is displayed. The target, slope,  $R^2$  value, Y-intercept, amplification efficiency, and error are displayed below the plot.

3. Confirm that the slope,  $R^2$  value, amplification efficiency, and error meet the analysis criteria.

For more information, see “Standard Curve Plot overview” on page 18.

4. Visually check that all unknown sample  $C_q$  values fall in the standard curve range.
5. In the **Well Table**, confirm that the  $C_q$  values of all replicate samples meet the analysis criteria.
6. In the **Replicate Group Table**, review the quantity mean and quantity SD if needed.

If the results do not meet the analysis criteria, troubleshoot using one of the following strategies:

- Omit wells, then reanalyze (see “Identify and omit outliers from standard curve analysis” on page 15).
- Repeat the plate run, adjusting plate file setup and analysis settings to improve results.

## Identify and omit outliers from standard curve 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 **Standard Curve** tab, in the plot pane, click on an outlier data point to highlight the well in the **Well Table**.  
Outlier wells can also be omitted in the **Quality Check** tab (see [? Help ▶ Help Contents](#)).
2. In the **Well Table**, select **Omit** in the row of the outlier well.
3. 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.

## (Optional) Edit standard curve analysis settings

Open the Standard Curve Analysis Module.

1. Click **Actions ▶ Standard Curve Analysis Setting**.
2. In the **General Setting** tab, edit the analysis settings if needed.

Standard Curve Analysis Option	Description
On Plate Standard Curves	Select to use the standard curve from the current data file. Click <b>Export</b> to export the standard curve.
External Standard Curves	Select to use a standard curve that was previously exported from another data file for analysis in the current data file. The two data files must be from the same instrument type, block type, and run method. <ul style="list-style-type: none"> <li>• Click <b>Import</b>, navigate to the standard curve file, then click <b>Open</b>.</li> <li>• Click <b>Delete</b> to delete an imported standard curve.</li> <li>• Click <b>Export</b> to export the standard curve.</li> </ul>

3. Click **Apply**.

The data is reanalyzed using the updated analysis settings.





# About standard curve analysis

## Overview of standard curve analysis

Standard curve analysis is used to determine absolute target quantity in samples.

For standard curve analysis, the software performs the following tasks.

1. The software measures amplification of the target in a standard dilution series and in test samples.
2. The software generates a standard curve using data from the standard dilution series.
3. The software uses the standard curve to interpolate the absolute quantity of the target in the test samples.

## Sample types for standard curve analysis

Standard curve analysis includes the following sample types for each target of interest. Each unique target requires its own standard curve.

Sample type (Type column in Samples table)	Sample description	Automatic target task assignment <sup>[1]</sup> (Task column in Targets table)
Standard	<p>A sample that contains known quantities or known relative quantities of the target</p> <ul style="list-style-type: none"><li>• For known quantities—quantify the target in the standard sample using an independent method</li><li>• For known relative quantities—generate a relative dilution series of the target standards</li></ul> <p><b>Note:</b> You must enter a quantity for each standard sample in the <b>Samples</b> table. Do not edit the quantity in the <b>Targets</b> table.</p>	Standard

Sample type (Type column in Samples table)	Sample description	Automatic target task assignment <sup>[1]</sup> (Task column in Targets table)
Unknown	Test sample	Unknown
Negative Control	Water or buffer No amplification of the target should occur in NTC wells.	Negative control

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

## Standard Curve Plot overview

The **Standard Curve Plot** displays the standard curve for samples designated as standards. The software calculates the quantity of a target in an unknown sample using the standard curve.

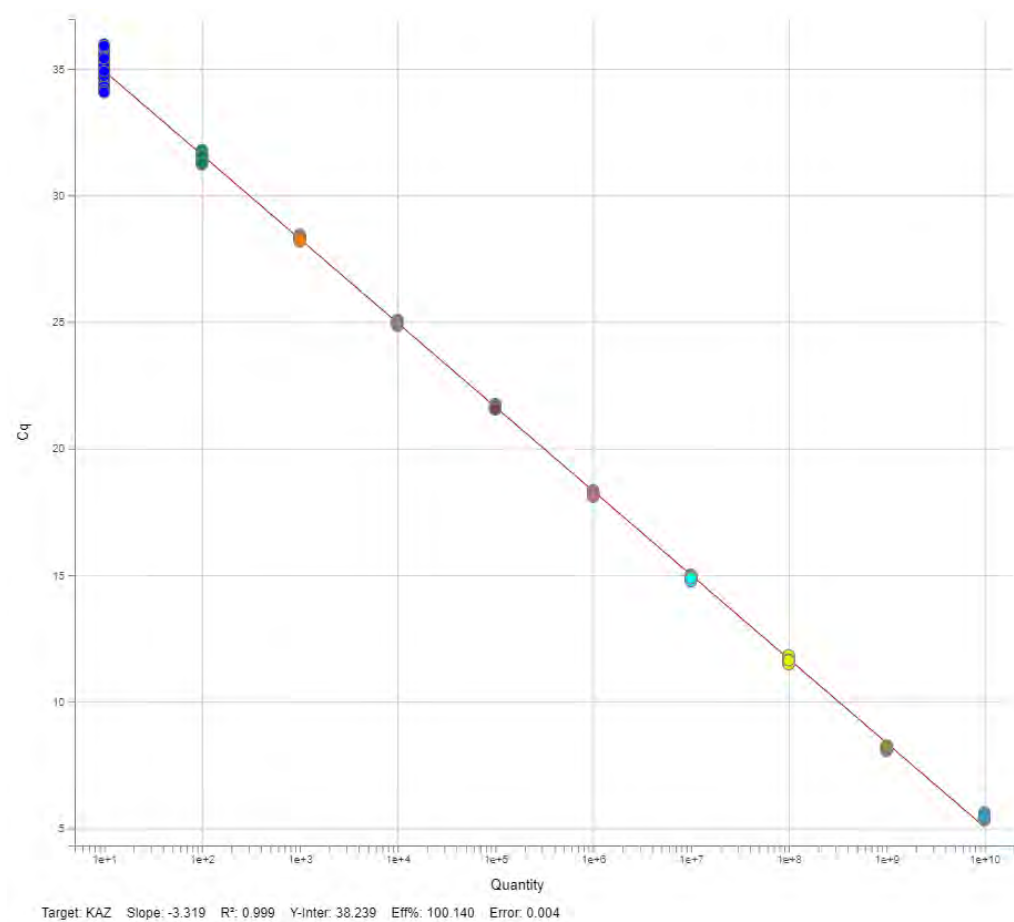


Figure 1 Example Standard Curve Plot

Table 1 Results or metrics to review in the Standard Curve Plot

Results or metrics	Description	Criteria for evaluation
Slope and amplification efficiency	The amplification efficiency is calculated using the slope of the regression line in the standard curve.	<p>A slope close to <math>-3.3</math> indicates optimal, 100% PCR amplification efficiency.</p> <p>Factors that affect amplification efficiency:</p> <ul style="list-style-type: none"> <li>• Improper design of the primer and probe</li> <li>• Range of standard quantities—For accurate and precise efficiency measurements, use a broad range of standard quantities, 5 to 6 logs (<math>10^5</math>- to <math>10^6</math>-fold).</li> <li>• Number of standard replicates—For accurate efficiency measurements, include replicates to decrease the effects of pipetting inaccuracies.</li> <li>• PCR inhibitors—PCR inhibitors and contamination in the reaction can reduce amplification efficiency.</li> <li>• Other possible factors: <ul style="list-style-type: none"> <li>– Component and properties of the reaction mix, such as salt content, DMSO, pH, etc.</li> <li>– Inaccurate sample or reagent pipetting</li> <li>– Improper analysis settings</li> <li>– Incorrect plate setup</li> </ul> </li> </ul>
$R^2$ value (correlation coefficient)	The $R^2$ value is a measure of the closeness of fit between the regression line and the individual $C_q$ data points of the standard reactions.	<ul style="list-style-type: none"> <li>• A value of 1.00 indicates a perfect fit between the regression line and the data points.</li> <li>• An <math>R^2</math> value <math>&gt; 0.99</math> is desirable.</li> </ul>
Error	<p>The standard error of the slope of the regression line in the standard curve.</p> <p>The error can be used to calculate a confidence interval (CI) for the slope and therefore the amplification efficiency.</p>	Acceptable value is determined by the analysis criteria.
$C_q$ values	$C_q$ is the PCR cycle number at which the fluorescence level meets the threshold.	<p>A <math>C_q</math> value <math>&gt; 8</math> and <math>&lt; 35</math> is desirable.</p> <ul style="list-style-type: none"> <li>• <math>C_q</math> value <math>&lt; 8</math>—There may be too much template in the reaction.</li> <li>• <math>C_q</math> value <math>&gt; 35</math>—There may be a low amount of target in the reaction; for <math>C_q</math> values <math>&gt; 35</math>, expect a higher standard deviation.</li> </ul>

# Documentation and support

## Related documentation

Document	Publication number
<i>QuantStudio™ Design and Analysis Software v2 User Guide</i>	MAN0018200
<i>QuantStudio™ 6 Pro and 7 Pro Real-Time PCR Systems User Guide</i>	MAN0018045
<i>QuantStudio™ 6 Pro and 7 Pro Real-Time PCR Systems Site Preparation Guide</i>	MAN0017992

## Customer and technical support

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

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

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

## Limited product warranty

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

## A

absolute quantity 17  
Amplification Plot, review 14  
analysis module, select 14

## C

compatible data file types 5  
Cq precision, improve 15

## L

limited product warranty 20

## M

MultiComponent Plot, review 15

## O

omit wells, Standard Curve 15

## P

passive reference, select 12  
perform analysis 14  
plate file  
    analysis type 8  
    create 8  
    save 12  
    set up 8, 9  
plate setup  
    samples 10  
    targets 10

## R

Raw Data Plot, review 16  
reagents, information 12  
related documentation 20  
review, Standard Curve Plot 18  
run method 9

## S

sample types, standard curve analysis 17  
samples 10  
standard curve  
    analysis 7  
    set up 11  
    workflow 7  
standard curve analysis  
    overview 17  
    sample types 17  
    samples 10  
    targets 10  
standard curve analysis module, about 5  
Standard Curve Plot  
    overview 18  
    review 14  
support, customer and technical 20

## T

target 10  
terms and conditions 20

## W

warranty 20

