## appliedbiosystems

## Custom Dye Calibration of Applied Biosystems<sup>™</sup> Real-Time PCR Instruments

Publication Number MAN0010765 Revision B.0

| About this guide   | 1  |
|--|----|
| Custom dyes overview   | 1  |
| Compatible instruments   | 2  |
| Workflow: Custom dye calibration                                       | 3  |
| Use a dilution series to determine an optimal custom dye concentration | 3  |
| Calibrate the custom dye   | 7  |
| Troubleshoot calibration failure                                       | 10 |
| Documentation and support  | 11 |

### About this guide

This document describes custom dye calibration procedures for:

- Interchangeable-block instruments (see "Compatible instruments" on page 2).
- Plate formats of 96-well Fast (0.1-mL), 96-well Standard (0.2-mL), and 384-well.

### **Custom dyes overview**

Applied Biosystems<sup>™</sup> real-time PCR instruments and systems can run assays designed with custom dyes. Custom dyes include:

- Dyes that are not manufactured by Thermo Fisher Scientific.
- Dyes or formulations of dyes that are not system dyes for the instrument.

To use a custom dye on the instrument, review the following requirements:

- Ensure that the custom dye excites and emits within the supported wavelength ranges for the instrument.
- Select a custom dye that does not overlap with other dyes used in the run.
- The custom dye must be attached to the 5' end of a short DNA oligonucleotide consisting of the first two bases of the probe sequence without a quencher at the 3' end.
- Calibrate the instrument for the custom dye.

Note: See the instrument maintenance guide for information about the following:

- · Supported wavelength ranges for the instrument
- · Descriptions of dyes (including their wavelengths) that are considered system dyes



# About dye<br/>calibrationThe software uses calibration data to characterize and distinguish the individual<br/>contribution of each dye in the total fluorescence signals collected by the instrument.

The software extracts a spectral profile for each dye standard, then produces a set of spectral profiles plotted as fluorescence versus filter.

The pass criteria for a dye calibration is that the dye spectra peak within the same filter as their group.

### **Compatible instruments**

The procedures in this guide are compatible with the following interchangeable-block real-time PCR instruments or systems. These procedures replace the custom dye calibration described in the user guides for the following instruments or systems:

| Instrument or system  | User guide  | Pub. No. |
|---|---|----------|
| ViiA <sup>™</sup> 7 Real-Time PCR System                            | Applied Biosystems <sup>™</sup> ViiA <sup>™</sup> 7 Real-Time PCR<br>System User Guide: Calibration, Maintenance,<br>Networking, and Security | 4442661  |
| QuantStudio <sup>™</sup> 6 Flex Real-Time PCR System                | QuantStudio <sup>™</sup> 6 and 7 Flex Real-Time PCR   | 4489821  |
| QuantStudio <sup>™</sup> 7 Flex Real-Time PCR System                | Systems Maintenance and Administration Guide  |          |
| QuantStudio <sup>™</sup> 12K Flex Real-Time PCR System              | QuantStudio <sup>™</sup> 12K Flex Real-Time PCR System<br>Maintenance and Administration Guide  | 4470689  |
| QuantStudio <sup>™</sup> Dx Real-Time PCR Instrument <sup>[1]</sup> | User guide or getting started guide for the<br>QuantStudio <sup>™</sup> Test Development Software   | _[2]     |

[1] In this document, for the QuantStudio<sup>™</sup> Dx Real-Time PCR Instrument, the term 'desktop software' refers to the QuantStudio<sup>™</sup> Test Development Software is for Research Use Only.

<sup>[2]</sup> See the user guide that corresponds to your version of QuantStudio<sup>™</sup> Test Development Software.

**Note:** Some of the listed compatible instruments can support TaqMan<sup>®</sup> Array Cards. For custom dye calibration procedures for array cards, contact your Thermo Fisher Scientific service representative.

### Workflow: Custom dye calibration

For each custom dye, determine the optimal dye concentration. Use this concentration to prepare all subsequent dye calibration plates.

Use a dilution series to determine an optimal custom dye concentration

Prepare a custom dye dilution plate (page 4)

Run the dilution plate as an experiment (page 5)

▼

Determine the optimal dye concentration (page 6)

Calibrate the custom dye using the optimal concentration

Create a custom dye calibration plate (page 7)

7

Add a new custom dye to the desktop software (page 7)

Perform a custom dye calibration (page 8)

# Use a dilution series to determine an optimal custom dye concentration

Custom dye dilution guidelines

Prepare a dilution series for each custom dye.

- Target several dye concentrations within a range of 100–2,000 nM.
- Choose a 2- or 3-fold difference in dilution points.
- Dispense an appropriate volume for the plate format.
  - 96-well Standard plate: 20 μL per well
  - 96-well Fast plate: 10 μL per well
  - 384-well plate: 10 µL per well
- Dilute the dye in buffer compatible with your master mix.
- (*Intercalating dyes only*) Add the appropriate amount of amplified PCR product to generate fluorescence.

# Prepare a custom dye dilution plate

**IMPORTANT!** Wear powder-free gloves throughout the procedure.

- 1. Prepare a 2- or 3-fold dilution series of the custom dye.
- **2.** Dispense aliquots of each dilution into the center of a reaction plate, then seal the plate.

A full plate is not needed. See the following figure for suggested replicates.





- **3.** Vortex the plate for 5 seconds, then centrifuge at 750 to  $1,000 \times g$  for 2 minutes.
- **4.** Confirm that the liquid in each well is at the bottom of the well and free of bubbles. If it is not, centrifuge the plate again.



**IMPORTANT!** Keep the bottom of the plate clean. Fluids and other contaminants on the bottom of the plate can contaminate the sample block and cause an abnormally high background signal.

### Run the dilution plate as an experiment

1. Load the plate into the instrument.

**IMPORTANT!** The instrument should be used by trained operators who have been warned of the moving parts hazard.

- 2. Set up a genotyping experiment in the desktop software.
  - a. In the **Home** tab, select **New Experiment Experiment Setup**.
  - **b.** In the **Properties** screen:
    - Select **Genotyping** as the experiment type.
    - At the bottom of the screen:
      - Deselect both Pre-PCR Read and Amplification.
      - Ensure that **Post-PCR Read** is selected (default).
    - (*Optional*) Enter or edit other experiment properties as needed.
  - **c.** In the **Define** and **Assign** screens, enter the dilution series information for the plate wells.
  - d. In the Run Method screen:
    - Enter the appropriate reaction volume.
    - Set the hold time to 2 minutes.
    - Set the hold temperature to the detection temperature used in your experiments.
       For TaqMan<sup>®</sup> Assays, we recommend 60°C.
    - Ensure the correct filter for the custom dye is selected. If you do not know the correct filter settings, select all the available filter combinations.



- **3.** Save the experiment, then start the run.
- 4. When the run ends, the EDS file automatically transfers to the desktop software.
- **5.** Unload the plate from the instrument.

**CAUTION!** PHYSICAL INJURY HAZARD. During instrument operation, the plate temperature can reach 100°C. Allow it to cool to room temperature before handling.

Note: If the instrument does not eject the tray arm, contact Support.

# Determine the optimal dye concentration

Review the dye signal data and select the dilution to use for dye calibrate.

- In the left menu pane, select Analysis > Raw Data Plot. This plot displays the raw fluorescence signal of each optical filter, for individual wells.
- **2.** For each replicate population of dilutions, select the wells in the plate layout to view in the plot.
- **3.** Examine the raw data to identify the wells yielding signals according to the ranges shown in the following table.

| Plate type | Acceptable signal range <sup>[1]</sup> |
|------------|--|
| 96-well    | 1,400,000 to 4,300,000                 |
| 384-well   | 400,000 to 1,200,000                   |

<sup>[1]</sup> Signal range for the optical filter where the dye is brightest.

- **4.** (*Optional*) Export the raw data, then calculate the average fluorescence value for each concentration.
- **5.** Select the lowest (optimal) dye concentration that falls within the acceptable signal range.

### Calibrate the custom dye

| Create  | а  | cus  | stom |
|---------|----|------|------|
| dye cal | ib | rati | on   |
| plate   |    |      |      |

**IMPORTANT!** Wear powder-free gloves while creating the dye plate.

Create a full plate of the custom dye diluted to the optimal concentration.

1. Dilute the custom dye to the optimal concentration in buffer.

Prepare an adequate volume, using an appropriate volume for the plate format.

- 96-well Standard plate: 20 μL per well
- 96-well Fast plate: 10 µL per well
- 384-well plate: 10 µL per well
- 2. Pipet the appropriate volume of the diluted custom dye to the plate wells.
- 3. Seal the plate.
- **4.** Vortex the plate for 5 seconds, then centrifuge at 750 to  $1,000 \times g$  for 2 minutes.
- **5.** Confirm that the liquid in each well is at the bottom of the well and free of bubbles. If it is not, centrifuge the plate again.



**IMPORTANT!** Keep the bottom of the plate clean. Fluids and other contaminants on the bottom of the plate can contaminate the sample block and cause an abnormally high background signal.

1. In the desktop software 🚰 Home tab, click 🧾 Instrument Console.

dye to the desktop 2. 1

Add a new custom

**2.** Select the instrument, then click **Manage Instrument**.

**Note:** If **Manage Instrument** is inactive, add your instrument (see the instrument maintenance guide).

- **3**. In the left menu pane, select **X Maintenance Dye**.
- 4. Select Custom Dye Calibration, then click Start Calibration.
- 5. Click New Dye, then in the Dye Library dialog box, click New.

6. Enter the custom dye information (see the following table), then click OK ► Close.

| Field/option | Action  |  |  |
|--------------|---|--|--|
| Name         | Enter a name for the custom dye.  |  |  |
|              | IMPORTANT!  |  |  |
|              | <ul> <li>Do not use a system dye name for a custom dye name.</li> <li>Dye names are spacing- and case-sensitive and cannot contain special characters.</li> </ul> |  |  |
| Wavelength   | Enter the wavelength at which the dye fluoresces.   |  |  |
| Туре         | Select:   |  |  |
|              | • <b>Reporter</b> —The dye works in conjunction with a quencher dye to report an increase of PCR product.   |  |  |
|              | • <b>Quencher</b> —The dye suppresses the fluorescence of a reporter dye until amplification of PCR product.  |  |  |
|              | • <b>Both</b> —The dye can be used as a reporter or quencher dye.   |  |  |

## Perform a custom dye calibration

### **1.** Load the plate into the instrument.

**IMPORTANT!** The instrument should be used by trained operators who have been warned of the moving parts hazard.

- 2. In the desktop software 🕂 Home tab, click 🧾 Instrument Console.
- 3. Select the instrument, then click Manage Instrument.

**Note:** If **Manage Instrument** is inactive, add your instrument (see the instrument maintenance guide).

- **4.** In the left menu pane, select **X Maintenance** > **Dye**.
- 5. Select Custom Dye Calibration, then click Start Calibration.
- 6. Select a custom dye from the dropdown list.
- 7. Enter the calibration temperature.

**Note:** The calibration temperature is the temperature at which the software will collect data. For TaqMan<sup>®</sup> reagents, we recommend using 60°C.

- 8. In the **Reagent Information** pane:
  - **a**. Enter the reagent information for the plate.
  - **b.** Select the checkbox to confirm that you loaded the plate and entered the reagent information.
  - c. Click Next.
- 9. Click Start Run.

**10.** When the run is complete, review the analysis status and dye spectra. Confirm that the dye signal peaked in the correct filter.

| Analysis status | Action   |
|-----------------|--|
| Passed          | Click Next.  |
| Caution         | See "Troubleshoot calibration failure" on page 10. |



Pass criteria—Signals from each well follow a uniform trend, and the dye peaks at the correct filter.

- 11. Click **Finish**, then follow the prompts to save the calibration results.
- **12.** Unload the plate from the instrument.



**CAUTION!** PHYSICAL INJURY HAZARD. During instrument operation, the plate temperature can reach 100°C. Allow it to cool to room temperature before handling.

Note: If the instrument does not eject the tray arm, contact Support.

## Troubleshoot calibration failure

| Observation   | Possible cause   | Recommended action  |
|---|--|---|
| Calibration result is <b>Caution</b>  | The plate was improperly prepared.   | <ul> <li>Ensure the following:</li> <li>The correct plate was used for the calibration performed.</li> <li>The plate was properly centrifuged.</li> <li>The plate was properly sealed.</li> <li>Consistent pipetting from well-to-well</li> </ul>   |
|   | The plate is damaged or contaminated.  | Check for damage, improper plate seal, or contamination.  |
| High fluorescence signal in<br>individual wells   | Signals that exceed the limit of<br>normal fluorescence may<br>indicate fluorescent<br>contaminants on the plate or<br>the sample block. | See the instrument maintenance guide for information about identifying contamination.   |
| High fluorescence signal<br>across the plate<br><i>(Custom dye calibrations only)</i>                                 | The dye concentration used to create the custom dye plate is too high.   | Create a new custom dye plate using the next<br>dye concentration lower than the optimal<br>concentration, then perform the calibration<br>again.   |
|   | A different temperature was<br>used for instrument calibration<br>than was used for determining<br>the optimal dye concentration.        | Ensure that the temperatures are consistent.  |
| Low fluorescence signal across<br>the plate<br><i>(Custom dye calibrations only)</i>                                  | A different temperature was<br>used for instrument calibration<br>than was used for determining<br>the optimal dye concentration.        | Ensure that the temperatures are consistent.  |
| Unexpected peaks or peaks<br>exceed maximum limit<br>• Spectra contain peaks that<br>are detected in more than        | Fluorescent contaminants on the sample block or on the plate.  | Perform a background calibration to detect<br>sample block contamination. If the background<br>calibration does not detect contamination, the<br>dye plate may be contaminated.   |
| <ul> <li>one filter.</li> <li>One or more raw spectra<br/>exceed the maximum limit<br/>for the instrument.</li> </ul> | <i>(Custom dye calibrations only)</i><br>Dye may be too concentrated.  | Prepare a new custom dye calibration plate but<br>decrease the concentration of the dye that<br>exceeds the detectable limit.   |
| One or more raw spectra are at<br>or below the detectable<br>threshold for the calibration                            | <ul> <li>The dye plate was<br/>centrifuged insufficiently.</li> <li>The dye plate contains old<br/>or insufficient reagents.</li> </ul>  | <ul> <li>Examine the wells of the plate.</li> <li>Liquid in the wells is not at the bottom of the wells—Centrifuge the plate for a longer time, then repeat the calibration.</li> <li>Liquid in the wells are equivalent in volume—The plate is improperly sealed and the reagents have evaporated. Discard the plate, then prepare and run a new plate.</li> <li>Examine the plate and the plate wells. If the plate appears to be normal, discard the plate,</li> </ul> |
|   |  | If the problem persists, contact Support.   |

| Observation  | Possible cause  | Recommended action  |
|--|---|---|
| One or more raw spectra are at<br>or below the detectable<br>threshold for the calibration | <i>(Custom dye calibrations only)</i><br>Dye may not be present at a<br>sufficient concentration. | Create a new custom dye plate using the next<br>dye concentration greater than the optimal<br>concentration, then perform the calibration<br>again. |
| Dye calibration passed, but the spectra did not peak in the correct filters                | The ROI calibration or<br>background calibration is<br>invalid.                                   | See the instrument maintenance guide for instructions on performing instrument calibrations.  |

### **Documentation and support**

| Customer and<br>technical support | <ul> <li>Visit thermofisher.com/support for the latest in services and support, including:</li> <li>Worldwide contact telephone numbers</li> <li>Product support, including: <ul> <li>Product FAQs</li> <li>Software, patches, and updates</li> <li>Training for many applications and instruments</li> </ul> </li> <li>Order and web support</li> <li>Product documentation, including: <ul> <li>User guides, manuals, and protocols</li> <li>Certificates of Analysis</li> <li>Safety Data Sheets (SDSs; also known as MSDSs)</li> <li>Note: For SDSs for reagents and chemicals from other manufacturers,</li> </ul> </li> </ul> |
|-----------------------------------|---|
| Limited product<br>warranty       | contact the manufacturer.<br>Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth<br>in the Life Technologies' General Terms and Conditions of Sale found on Life<br>Technologies' website at www.thermofisher.com/us/en/home/global/<br>terms-and-conditions.html. If you have any questions, please contact Life<br>Technologies at www.thermofisher.com/support.  |

### Symbols that may be displayed on the instrument, in the software, or in this guide

| Symbol | Description                  | Symbol | Description                                |
|--------|------------------------------|--------|--|
|        | MANUFACTURER                 | ~      | DATE OF MANUFACTURE                        |
| REF    | CATALOG NUMBER               | SN     | SERIAL NUMBER                              |
| Ĩ      | CONSULT INSTRUCTIONS FOR USE | Â      | CAUTION, CONSULT ACCOMPANYING<br>DOCUMENTS |



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

#### 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. MAN001765

| Revision | Date           | Description  |
|----------|----------------|--|
| B.0      | 18 April 2018  | <ul> <li>Update branding, general formatting, and images.</li> <li>Streamline content and phrasing.</li> <li>Clasify custom dvo consolution and extract setup for determining entimal system dvo concentration.</li> </ul> |
|          |                | <ul> <li>Add applicable instruments, troubleshooting items, and recommended volumes for plate formats.</li> </ul>  |
| A.0      | 27 August 2014 | New document.  |

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

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

©2018 Thermo Fisher Scientific Inc. All rights reserved.

#### thermofisher.com/support | thermofisher.com/askaquestion



thermofisher.com