

# User Bulletin

## Applied Biosystems StepOne™ and StepOnePlus™ Real-Time PCR Systems

March 2008

**SUBJECT: System Software Patch Provides Support for Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub>™ 1-Step Kits**

### In This User Bulletin

This user bulletin describes how to:

[Download and Install the StepOne™ Software v2.0.2 Patch . . . . . 3](#)

[Use Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub>™ 1-Step Kits . . . . . 7](#)

### About the Software Patches

The StepOne™ Software v2.0.2 patch:

- Allows you to use the Design Wizard, Advanced Setup, and QuickStart features to design and set up quantitation experiments using the Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub>™ 1-Step Kits. With the patch installed, the StepOne™ software uses the appropriate default run methods for the Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub>™ 1-Step Kits.
- Displays the relevant reagents in the Materials List of the StepOne™ software when you select SYBR® Green reagents or TaqMan® reagents and the applicable Fast or Standard ramp speed.

### Before You Begin

The following table lists the versions of software that must be present on your StepOne™ or StepOnePlus™ system computer before you can install the update. Before installing the update, make sure that your computer is running the appropriate version of the software.

System	The computer must have installed...
StepOne™ system	<ul style="list-style-type: none"><li>• StepOne™ Software v2.0</li><li>or</li><li>• StepOne™ Software v2.0.1</li></ul>
StepOnePlus™ system	

## Related Products

### Table 1 Fast SYBR<sup>®</sup> Green Master Mix Contents

Kit/Contents	20-μL Rxns/Kit	Part Number
Fast SYBR <sup>®</sup> Green Master Mix (2X), 1 mL	100	4385610
Fast SYBR <sup>®</sup> Green Master Mix (2X), 5 mL	500	4385612
Fast SYBR <sup>®</sup> Green Master Mix (2X), 2 × 5 mL	1000	4385616
Fast SYBR <sup>®</sup> Green Master Mix (2X), 5 × 5 mL	2500	4385617
Fast SYBR <sup>®</sup> Green Master Mix (2X), 10 × 5 mL	5000	4385618
Fast SYBR <sup>®</sup> Green Master Mix (2X), 50 mL	5000	4385614

### Table 2 Power SYBR<sup>®</sup> Green RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Contents

Kit/Contents	50-μL Rxns/Kit	Part Number
<ul style="list-style-type: none"> <li>Power SYBR<sup>®</sup> Green RT-PCR Mix (2X), 1 mL</li> <li>RT Enzyme Mix (125X), 20 μL</li> </ul>	40	4391178
<ul style="list-style-type: none"> <li>Power SYBR<sup>®</sup> Green RT-PCR Mix (2X), 5 mL</li> <li>RT Enzyme Mix (125X), 80 μL</li> </ul>	200	4389986

### Table 3 TaqMan<sup>®</sup> RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Contents

Kit/Contents	50-μL Rxns/Kit	Part Number
<ul style="list-style-type: none"> <li>TaqMan<sup>®</sup> RT-PCR Mix (2X), 1 mL</li> <li>TaqMan<sup>®</sup> RT Enzyme Mix (40X), 50 μL</li> </ul>	40	4392653
<ul style="list-style-type: none"> <li>TaqMan<sup>®</sup> RT-PCR Mix (2X), 5 mL</li> <li>TaqMan<sup>®</sup> RT Enzyme Mix (40X), 250 μL</li> </ul>	200	4382938
<ul style="list-style-type: none"> <li>TaqMan<sup>®</sup> RT-PCR Mix (2X), 10 × 5 mL</li> <li>TaqMan<sup>®</sup> RT Enzyme Mix (40X), 10 × 250 μL</li> </ul>	2000	4392656

## Related Documentation

- Access the StepOne<sup>™</sup> Software Help by clicking [?](#) or pressing **F1**.
- Applied Biosystems StepOne<sup>™</sup> and StepOnePlus<sup>™</sup> Real-Time PCR Systems Getting Started Guide for Standard Curve Experiments* (PN 4376784)
- Applied Biosystems StepOne<sup>™</sup> and StepOnePlus<sup>™</sup> Real-Time PCR Systems Getting Started Guide for Relative Standard Curve and Comparative C<sub>T</sub> Experiments* (PN 4376785)
- Fast SYBR<sup>®</sup> Green Master Mix Protocol* (PN 4385372)
- Fast SYBR<sup>®</sup> Green Master Mix Quick Reference Card* (PN 4385371)
- Power SYBR<sup>®</sup> Green RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Protocol* (PN 4391003)
- Power SYBR<sup>®</sup> Green RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Quick Reference Card* (PN 4391112)
- TaqMan<sup>®</sup> RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Protocol* (PN 4393463)
- TaqMan<sup>®</sup> RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit Quick Reference Card* (PN 4393464)

# Download and Install the StepOne™ Software v2.0.2 Patch

Perform the following procedure to download and install the StepOne™ Software v2.0.2 patch for Applied Biosystems StepOne™ and StepOnePlus™ Fast Real-Time PCR Systems.

## Download the Patch

1. Go to the StepOne™ or StepOnePlus™ system website:
  - <http://info.appliedbiosystems.com/stepone>
  - <http://info.appliedbiosystems.com/steponeplus>
2. In the Software Download page, click **Download StepOne Software v2.0.2 Patch**.

### Applied Biosystems introduces the 96-well StepOnePlus™ Real-Time PCR System

#### 48-Well StepOne™ System

- Surprisingly affordable
- 48 wells for lower throughput
- 3 colors for basic applications research
- Easily upgradeable to the StepOnePlus™ System

#### New! 96-Well StepOnePlus™ System

- 96-wells for higher throughput
- 4 colors for more versatility
- Same easy to use software as the StepOne™ System

**The Affordable 48-Well StepOne™ System Offers a Seamless Upgrade to the 96-Well StepOnePlus™ System with Easy-to-Use Software...A Step Above the Rest.**  
 The StepOne and StepOnePlus Real-Time PCR Systems deliver remarkable plug-and-play convenience and uncompromising performance. Both systems support the full range of TaqMan® assays while providing ease of use and a small footprint. Each system offers intuitive software, fast runs (<40 minutes) and standard runs (<2 hours), and are configured for PC-free or networked operation. The factory calibrated StepOne and StepOnePlus Systems make it easy to step up to high performance real-time PCR.

[Download Brochure](#)
[View Software Demo](#)
[Request more Information](#)

#### Software Demo

- [Software Demo](#)

---

#### Product Information

- [StepOne/StepOnePlus System Brochure](#)
- [StepOne/StepOnePlus System Specification Sheet](#)
- [StepOne Product Page](#)
- [StepOnePlus Product Page](#)
- [Computer Specifications](#)

---

#### Software Download

- [Download StepOne Software v2.0](#)
- [Download StepOne Software v2.0.2 Patch](#)

2


3. In the Software Downloads page, click the icon (**PC**) for the **StepOne Software v2.0.2 Patch for StepOne and StepOnePlus Systems**.

## Software Downloads


StepOne and StepOnePlus Real-Time PCR Systems

Title	Software	Size	Related Documentation	PDF Size
StepOne Software v2.0.2 Patch for StepOne and StepOnePlus Systems	<b>PC</b>	25.4 MB	Read Me File	200k
Power SYBR Green RNA-to-Ct 1-Step 48-Well Template for StepOne Systems	<b>PC</b>	1k	Read Me File	121k
Power SYBR Green RNA-to-Ct 1-Step 96-Well Template for StepOnePlus Systems	<b>PC</b>	1k	Read Me File	121k
TaqMan RNA-to-Ct 48-Well 1-Step Template for StepOne Systems	<b>PC</b>	1k	Read Me File	121k
TaqMan RNA-to-Ct 96-Well 1-Step Template for StepOnePlus Systems	<b>PC</b>	1k	Read Me File	121k
Fast SYBR Green 48-Well Template for StepOne Systems	<b>PC</b>	1k	Read Me File	121k

3















4. In the File Download dialog box, click **Save**.
5. In the Save As dialog box, select a folder in which to save the patch, click **Save**, then click **Close**.
6. In the Software Downloads page, download the Read Me file for the patch:
  - a. Click the PDF () icon for the Read Me file.
  - b. In the File Download dialog box, click **Save**.
  - c. Select a folder in which to save the file, click **Save**, then click **Close**.
7. Continue to [“Download the Template”](#) below.

## Download the Template


1. In the Software Downloads page, click the appropriate icon () for your system:
  - For StepOne<sup>™</sup> systems, select one of the following:
    - **Fast SYBR Green 48-Well Template**
    - **Power SYBR Green RNA-to-C<sub>T</sub> 1-Step 48-Well Template**
    - **TaqMan RNA-to-C<sub>T</sub> 48-Well 1-Step Template**
  - For StepOnePlus<sup>™</sup> systems, select one of the following:
    - **Fast SYBR Green 96-Well Template**
    - **Power SYBR Green RNA-to-C<sub>T</sub> 96-Well 1-Step Template**
    - **TaqMan RNA-to-C<sub>T</sub> 96-Well 1-Step Template**

## Software Downloads

StepOne and StepOnePlus Real-Time PCR Systems

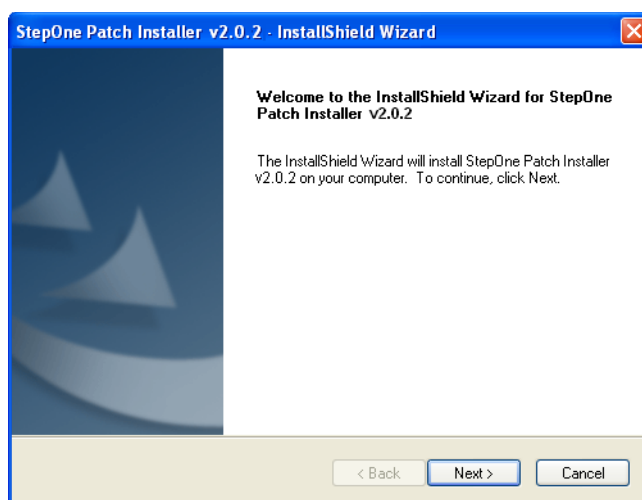
Title	Software	Size	Related Documentation	PDF Size
StepOne Software v2.0.2 Patch for StepOne and StepOnePlus Systems		25,4 MB	Read Me File	 208k
Power SYBR Green RNA-to-C <sub>T</sub> 1-Step 48-Well Template for StepOne Systems		1k	Read Me File	 121k
Power SYBR Green RNA-to-C <sub>T</sub> 1-Step 96-Well Template for StepOnePlus Systems		1k	Read Me File	 121k
TaqMan RNA-to-C <sub>T</sub> 48-Well 1-Step Template for StepOne Systems		1k	Read Me File	 121k
TaqMan RNA-to-C <sub>T</sub> 96-Well 1-Step Template for StepOnePlus Systems		1k	Read Me File	 121k
Fast SYBR Green 48-Well Template for StepOne Systems		1k	Read Me File	 121k
Fast SYBR Green 96-Well Template for StepOnePlus Systems		1k	Read Me File	 121k

1

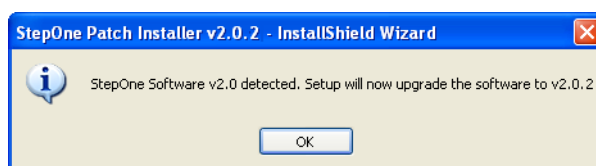
2. In the File Download dialog box, click **Save**.
3. In the Save As dialog box, select a folder in which to save the template zip file, click **Save**, then click **Close**.
4. Unzip the file.
5. In the Software Downloads page, download the corresponding Read Me file for the template:
  - a. Click the PDF icon () for the Read Me file.
  - b. In the File Download dialog box, click **Save**.
  - c. Select a folder in which to save the Read Me file, click **Save**, then click **Close**.
6. Continue to [“Install the Patch”](#) on page 5.

## Install the Patch

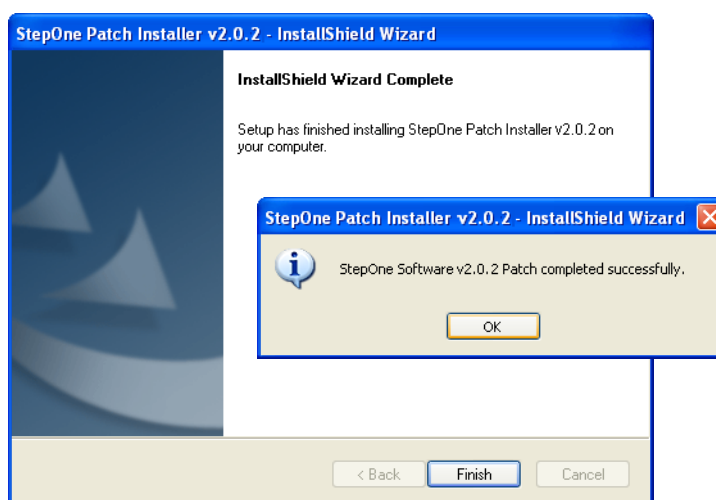
1. Double-click the **StepOne\_Software\_v2.0.2\_Patch.exe** file that you downloaded to your computer.
2. In the Welcome page of the InstallShield Wizard, click **Next**.



3. In the License Agreement page, click **Yes** to accept the agreement.  
**IMPORTANT!** You must accept the license agreement to install the patch.
4. After the installer detects StepOne™ Software v2.0 or v2.0.1, click **OK**.



5. After the successful installation, click **OK**, then click **Finish**.






6. Continue to [“Install the Template” on page 6](#).

## Install the Template

1. Transfer the template file from your computer to a USB drive:
  - a. Connect the USB drive to one of the USB ports on your computer.
  - b. Copy the **SYBR\_GREEN\_FST.ed**s, **SYBR\_RNA\_TO\_CT.ed**s, or **TQMN\_RNA\_TO\_CT.ed**s file from the computer to the USB drive.
2. Remove the USB drive from your computer, then connect the drive to the USB port of your StepOne™ or StepOnePlus™ instrument.



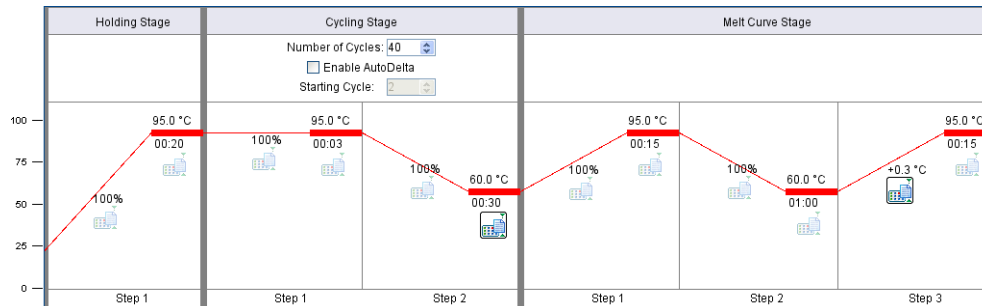
3. Touch the StepOne™ or StepOnePlus™ instrument touchscreen to awaken it.  
**Note:** If the touchscreen does not display the Main Menu screen, touch . If the instrument is powered off, touch .
4. Wait for the USB icon to appear on the touchscreen.  
**Note:** If your StepOne™ or StepOnePlus™ instrument cannot find the USB drive, remove the USB drive, then repeat [steps 2 through 4](#). If the instrument fails again to find the USB drive, repeat [steps 1 through 4](#) using another USB drive.
5. Transfer the template to the instrument:
  - a. In the Main Menu, touch **Browse/New Experiments**.
  - b. Touch  **Folders**, then touch **USB**.
  - c. In the Browse screen, touch **SYBR\_GREEN\_FST.ed**s, **SYBR\_RNA\_TO\_CT.ed**s, or **TQMN\_RNA\_TO\_CT.ed**s, then touch **Copy**.
  - d. In the Save Experiment screen, navigate to the Template folder, then touch **Save & Exit**.
6. When prompted that the data were transferred successfully, touch **OK**.
7. Remove the USB drive from the instrument.

# Use Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kits

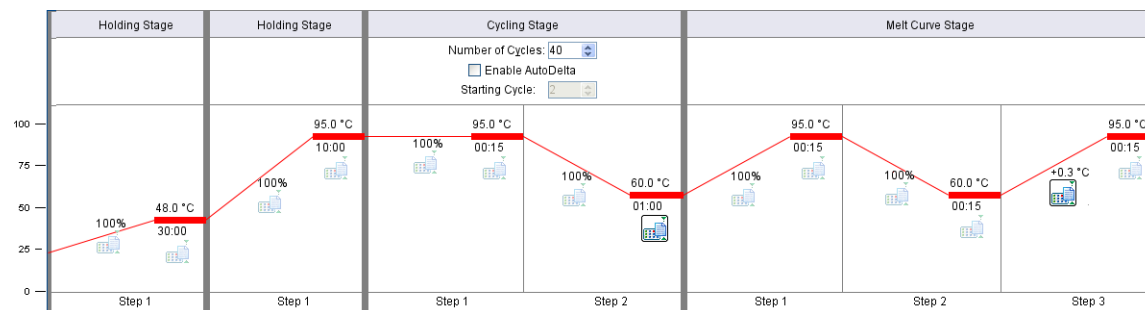
## StepOne<sup>™</sup> and StepOnePlus<sup>™</sup> System Thermal Profiles

After you install the updates, you can use the StepOne<sup>™</sup> Software v2.0.2 or the instrument touchscreen to perform quantitation experiments using the Fast SYBR® Green Master Mix and RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kits. When performing a run, make sure that the run method displays the thermal profile appropriate for your chemistry.

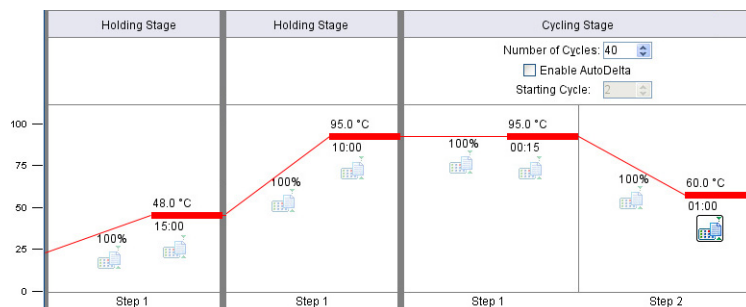
### Fast SYBR® Green Master Mix



### Power SYBR® Green RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit



### TaqMan® RNA-to-C<sub>T</sub><sup>™</sup> 1-Step Kit



© Copyright 2008, Applied Biosystems. All rights reserved.

**For Research Use Only. Not for use in diagnostic procedures.**

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

APPLIED BIOSYSTEMS DISCLAIMS ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL APPLIED BIOSYSTEMS BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

**TRADEMARKS:**

Applera, Applied Biosystems, and AB (Design) are registered trademarks, and RNA-to-C<sub>T</sub>, StepOne, and StepOnePlus are trademarks of Applera Corporation or its subsidiaries in the U.S. and/or certain other countries.

SYBR is a registered trademark of Molecular Probes, Inc.

TaqMan is a registered trademark of Roche Molecular Systems, Inc.

All other trademarks are the sole property of their respective owners.

**Worldwide Sales and Support**

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at [www.appliedbiosystems.com](http://www.appliedbiosystems.com).

Applera is committed to providing the world's leading technology and information for life scientists. Applera Corporation consists of the Applied Biosystems and Celera Genomics businesses.

**Headquarters**

850 Lincoln Centre Drive  
Foster City, CA 94404 USA  
Phone: +1 650.638.5800  
Toll Free (In North America): +1 800.345.5224  
Fax: +1 650.638.5884