

User Bulletin

Applied Biosystems RapidFinder™ Express Software v1.0
and SDS Software v1.4.1

January 2009

SUBJECT: Installing Applied Biosystems RapidFinder™ Express Software v1.0 and SDS Software v1.4.1

Overview The steps for installing a new Applied Biosystems 7500 Fast system with RapidFinder Express Software v1.0 and SDS Software v1.4.1 are provided in the *Applied Biosystems 7300/7500/7500 Fast Real-Time PCR System Installation and Maintenance Guide* (PN 4378657):

1. Plan the installation.
2. Uncrate and place the instrument.
3. Check the shipped materials. MicroSEQ® pathogen detection kits are not provided with the system when you purchase the RapidFinder™ Express Software.
4. Install the software using the Real-Time System Setup Wizard CD.
5. Perform a 7500 Fast Real-Time PCR System (ROI) calibration.
6. Perform a background calibration.
7. Perform an optical calibration.
8. Perform a pure-dye calibration.
9. Perform an RNase P verification run.

In This User Bulletin This user bulletin addresses requirements that you should meet before you install the software using the Real-Time System Setup Wizard CD (Step 4 above).

Minimum computer requirements

The minimum computer requirements for a 7500 Fast System with RapidFinder Express Software v1.0 and SDS Software v1.4.1 are as follows:

- 1 gigahertz
- 256 megabytes RAM
- Pentium III
- 20GB hard drive
- Microsoft Windows XP professional SP2

Computer resolution requirements

Set the Screen Resolution to a minimum of 1024 by 768 pixels with True Color (32 bit) (Control Panel ▶ Display ▶ Settings).

To install a new 7500 Fast system

RapidFinder Express Software v1.0 and SDS Software v1.4.1 are automatically installed on a new system during 7500 Fast installation. Insert the Real-Time System Setup Wizard CD into the computer CD drive and follow the instructions in the Real-Time System Setup Wizard.

To upgrade an existing 7500 Fast system

To upgrade your system:

1. Navigate to Applied Biosystems ▶ SDS Documents. Archive your SDS data.
2. On your control panel, click **Add/Remove Programs**. Uninstall the current version of SDS software.
3. Insert the Real-Time System Setup Wizard CD into the computer CD drive.
4. Select the **Install the SDS Software** option, then follow the instructions in the Real-Time Systems Setup Wizard.

© Copyright 2009, 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. This document is believed to be complete and accurate at the time of publication. In no event shall Applied Biosystems be liable for incidental, special, multiple, or consequential damages in connection with or arising from the use of this document.

NOTICE TO PURCHASER: LICENSE DISCLAIMER

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 Applied Biosystems, either expressly or by estoppel.

TRADEMARKS:

Applied Biosystems, AB (Design), and MicroSEQ are registered trademarks and RapidFinder is a trademark of Applied Biosystems or its subsidiaries in the US and/or certain other countries.

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.

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

1/2009