

Applied Biosystems ViiA™ 7 Real-Time PCR System Site Preparation Guide

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

Information in this document is subject to change without notice.

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. TO THE FULLEST EXTENT ALLOWED BY LAW, 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, WHETHER OR NOT FORESEEABLE AND WHETHER OR NOT APPLIED BIOSYSTEMS IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

Dell is a trademark of Dell Inc.

Milli-Q is a trademark of Millipore Corporation.

Twister is a trademark of Caliper Life Sciences Inc.

© 2010 Life Technologies Corporation. All rights reserved.

Part Number 4445302 Rev. B
12/2010

Contents

| | | |
|------------------|---|-----------|
| | About This Guide | 7 |
| | About this document | 7 |
| | Purpose of this guide | 7 |
| | Audience | 7 |
| | User attention words | 7 |
| | Safety conventions used in this document | 8 |
| | Safety alert words | 8 |
| | Definitions | 8 |
| | Examples | 8 |
| | Safety information | 10 |
| | Safety labels on instruments | 10 |
| | How to obtain support | 11 |
| | How to obtain more information | 12 |
| | Related documentation | 12 |
| | Send us your comments | 12 |
| CHAPTER 1 | Site Preparation Tasks | 13 |
| | Instrument overview | 14 |
| | Installation overview | 14 |
| | Site preparation schedule | 14 |
| | Site preparation process | 15 |
| | Assign personnel | 16 |
| | Laboratory safety representative | 16 |
| | Tasks and personnel | 16 |
| | Select the site | 18 |
| | Space requirements | 19 |
| | System components | 19 |
| | Setup requirements | 19 |
| | Layout requirements | 20 |
| | Dimensions and weights | 21 |
| | Required clearances | 21 |
| | Environmental requirements | 22 |
| | Altitude | 22 |
| | Temperature and humidity requirements | 22 |
| | Pollution | 22 |
| | Environmental conditions for storage and transportation | 22 |

| | |
|--|---------------|
| Instrument hot-air exhaust venting | 22 |
| Waste collection requirements | 23 |
| Collecting liquid | 23 |
| Electrical requirements | 23 |
| System electrical requirements | 23 |
| Electrical protective devices | 24 |
| Computer requirements | 25 |
| Antivirus software requirements | 25 |
| Network requirements | 25 |
| Safety and materials | 26 |
| Safety practices | 26 |
| Required safety equipment | 26 |
| Materials for installation | 27 |
| Materials for routine operation | 27 |
| Receiving and inspecting the system | 28 |
| Shipped contents | 28 |
| Shipping list | 28 |
| Inspect shipping containers for damage | 28 |
| Unpack and store the Installation Kit | 28 |
| Inspecting shipping containers for damage | 29 |
| Moving the crated instrument to the laboratory | 29 |
| Moving schedule | 29 |
| During installation | 30 |
| Operator training | 30 |
| CHAPTER 2 Checklists | 31 |
| Overview | 32 |
| Personnel checklist | 32 |
| Space and layout checklist | 32 |
| Environmental checklist | 33 |
| Electrical checklist | 33 |
| Computer checklist | 33 |
| Safety checklist | 34 |
| Materials checklist | 34 |
| System receipt and inspection checklist | 35 |
| Moving the crated instrument checklist | 35 |
| APPENDIX A Safety and EMC Compliance Information | 37 |
| Symbols on instruments | 38 |
| Electrical symbols on instruments | 38 |
| Safety symbols | 38 |
| Environmental symbols on instruments | 39 |

| | |
|--|-----------|
| Location of safety symbols | 40 |
| General instrument safety | 41 |
| Moving and lifting the instrument | 41 |
| Moving and lifting stand-alone computers and monitors | 41 |
| Operating the instrument | 41 |
| Cleaning or decontaminating the instrument | 42 |
| Chemical safety | 42 |
| Chemical hazard warning | 42 |
| About SDSs | 42 |
| Obtain an SDS | 42 |
| Chemical safety guidelines | 43 |
| Chemical waste safety | 44 |
| Chemical waste hazard | 44 |
| Chemical waste safety guidelines | 44 |
| Waste disposal | 45 |
| Electrical safety | 45 |
| Power | 45 |
| Overvoltage rating | 45 |
| Fuses | 46 |
| Bar code scanner laser safety | 46 |
| Physical hazard safety | 46 |
| Ultraviolet light | 46 |
| Moving parts | 46 |
| Biological hazard safety | 47 |
| General biohazard | 47 |
| Workstation safety | 47 |
| Safety and electromagnetic compatibility (EMC) standards | 48 |
| U.S. and Canadian safety standards | 48 |
| Canadian EMC standard | 48 |
| European safety and EMC standards | 48 |
| Australia and New Zealand EMC standards | 49 |
| Index | 51 |

About This Guide

About this document

Purpose of this guide

The *Applied Biosystems ViiA™ 7 Real-Time PCR System Site Preparation Guide* provides the information you need to fully prepare your site for the arrival and installation of the system. Complete preparation helps ensure a smooth installation process, as well as correct and safe instrument operation.

Audience

This guide is intended for the personnel who will schedule, manage, and perform the tasks required to prepare your site for installation of the *Applied Biosystems ViiA™ 7 Real-Time PCR System*.

User attention words

Two user attention words appear in Applied Biosystems user documentation. Each word implies a particular level of observation or action as described below:

Note: – Provides information that may be of interest or help but is not critical to the use of the product.

IMPORTANT! – Provides information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.

Examples of the user attention words appear below:

Note: For additional documentation, see [“How to obtain support”](#) on [page 11](#).

IMPORTANT! To verify your client connection to the database, you need a valid user ID and password.

Safety conventions used in this document

Safety alert words

Four safety alert words appear in Applied Biosystems user documentation at points in the document where you need to be aware of relevant hazards. Each alert word — **IMPORTANT**, **CAUTION**, **WARNING**, **DANGER** — implies a particular level of observation or action, as defined below.

Definitions

IMPORTANT! – Indicates information that is necessary for proper instrument operation, accurate chemistry kit use, or safe use of a chemical.



CAUTION! – Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.



WARNING! – Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.



DANGER! – Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury. This signal word is to be limited to the most extreme situations.

Except for IMPORTANTs, each safety alert word in an Applied Biosystems document appears with an open triangle figure that contains a hazard symbol. *These hazard symbols are identical to the hazard symbols that are affixed to Applied Biosystems instruments (see “Safety symbols” on page 38).*

Examples

The following examples show the use of safety alert words:

IMPORTANT! You must create a separate sample entry spreadsheet for each 384-well plate.



CAUTION! Do not install on the computer additional software other than antivirus software. Changes to the configured software could void the instrument warranty and cause the system to be nonoperational.



WARNING! Physical Injury Hazard. Using the instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument







DANGER! ELECTRICAL HAZARD. Failure to ground the instrument properly can lead to an electrical shock. Ground the instrument according to the provided instructions.

Safety information

Safety labels on instruments

The following CAUTION, WARNING, and DANGER statements may be displayed on Applied Biosystems instruments in combination with the safety symbols described in the preceding section.

| Hazard symbol | English | Français |
|---|---|---|
|  | CAUTION! Hazardous chemicals. Read the Safety Data Sheets (SDSs) before handling. | ATTENTION! Produits chimiques dangereux. Lire les fiches techniques de sûreté de matériels avant toute manipulation de produits. |
| | CAUTION! Hazardous waste. Refer to SDS(s) and local regulations for handling and disposal. | ATTENTION! Déchets dangereux. Lire les fiches techniques de sûreté de matériels et la réglementation locale associées à la manipulation et l'élimination des déchets. |
|  | WARNING! Hot lamp. | AVERTISSEMENT! Lampe brûlante. |
| | WARNING! Hot. Do not remove lamp until 15 min after disconnecting supply. | AVERTISSEMENT! Lampe brûlante, après avoir déconnecté le câble d'alimentation de l'appareil, attendre environ 15 minutes avant d'effectuer un remplacement de la lampe. |
| | WARNING! Hot. Replace lamp with an Applied Biosystems lamp. | AVERTISSEMENT! Composants brûlants. Remplacer la lampe par une lampe Applied Biosystems. |
| | CAUTION! Hot surface. | ATTENTION! Surface brûlante. |
|  | DANGER! High voltage. | DANGER! Haute tension. |
| | WARNING! To reduce the chance of electrical shock, do not remove covers that require tool access. No user-serviceable parts are inside. Refer servicing to Applied Biosystems qualified service personnel. | AVERTISSEMENT! Pour éviter les risques d'électrocution, ne pas retirer les capots dont l'ouverture nécessite l'utilisation d'outils. L'instrument ne contient aucune pièce réparable par l'utilisateur. Toute intervention doit être effectuée par le personnel de service qualifié venant de chez Applied Biosystems. |
|  | CAUTION! Moving parts. Crush/pinch hazard. | ATTENTION! Pièces en mouvement, risque de pincement et/ou d'écrasement. |

How to obtain support

For the latest services and support information for all locations, go to www.appliedbiosystems.com, then click the link for **Support**.

At the Support page, you can:

- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support
- Order Applied Biosystems user documents, certificates of analysis, and other related documents
- Order Safety Data Sheets (SDSs) for any chemicals supplied by Applied Biosystems or Ambion (available to you free 24 hours a day).
- Download PDF documents
- Obtain information about customer training
- Download software updates and patches

In addition, the Support page provides access to worldwide telephone and fax numbers to contact Applied Biosystems Technical Support and Sales facilities.

How to obtain more information

Related documentation

The following related documents are part of your shipment:

- **Applied Biosystems ViiA™ 7 Real-Time PCR System User Guide** – The user guide provides step-by-step instructions for preparing and analyzing a sample. It is designed to help you learn how to use the system.
- **Applied Biosystems ViiA™ 7 Real-Time PCR System Quick Reference Card** – The Quick Reference Card contains a flowchart on how to run the samples and system, a table of maintenance tasks, and a software reference guide.
- **Applied Biosystems ViiA™ 7 Real-Time PCR System Getting Started Guide** – The Getting Started Guide is a tutorial for performing your experiments.

Portable document format (PDF) versions of the User Guide and Quick Reference Card are also available on the Software CD.

Note: To view the user documentation, you can use the Adobe® Acrobat® Reader® software available from www.adobe.com.

Note: For additional documentation, see “How to obtain support” on page 11.

Send us your comments

Applied Biosystems welcomes your comments and suggestions for improving its user documents. You can e-mail your comments to:

techpubs@appliedbiosystems.com

IMPORTANT! The e-mail address above is only for submitting comments and suggestions relating to documentation. To order documents, download PDF files, or for help with a technical question, go to www.appliedbiosystems.com, then click the link for **Support**. (See “How to obtain support”).

1

Site Preparation Tasks

| | |
|--|----|
| ■ Instrument overview | 14 |
| ■ Installation overview | 14 |
| ■ Assign personnel | 16 |
| ■ Select the site | 18 |
| ■ Space requirements | 19 |
| ■ Environmental requirements | 22 |
| ■ Waste collection requirements | 23 |
| ■ Electrical requirements | 23 |
| ■ Computer requirements | 25 |
| ■ Safety and materials | 26 |
| ■ Materials for routine operation | 27 |
| ■ Receiving and inspecting the system | 28 |
| ■ Moving the crated instrument to the laboratory | 29 |
| ■ During installation | 30 |

Instrument overview

The instrument uses fluorescent-based polymerase chain reaction (PCR) reagents to provide:

- Quantitative detection of target nucleic acid sequences (targets) using real-time analysis.
- Qualitative detection of targets using post-PCR (endpoint) analysis.
- Qualitative analysis of the PCR product (achieved by melt-curve analysis that occurs post-PCR).

The instrument is to be used only by technologists trained in laboratory techniques and procedures and in the use of the instrument. The customer is responsible for validation of any assays, and compliance with any regulatory requirements that pertain to their procedures and uses of the instrument.

Installation overview

Before an Applied Biosystems service representative arrives to install the instrument and associated devices, prepare your site for the installation according to the instructions in this chapter. For checklists, see [“Checklists” on page 31](#).

IMPORTANT! If site preparation tasks are not complete when the Applied Biosystems service representative arrives, the scheduled installation may be postponed.

Site preparation schedule

To minimize the time between the shipment arrival and installation:

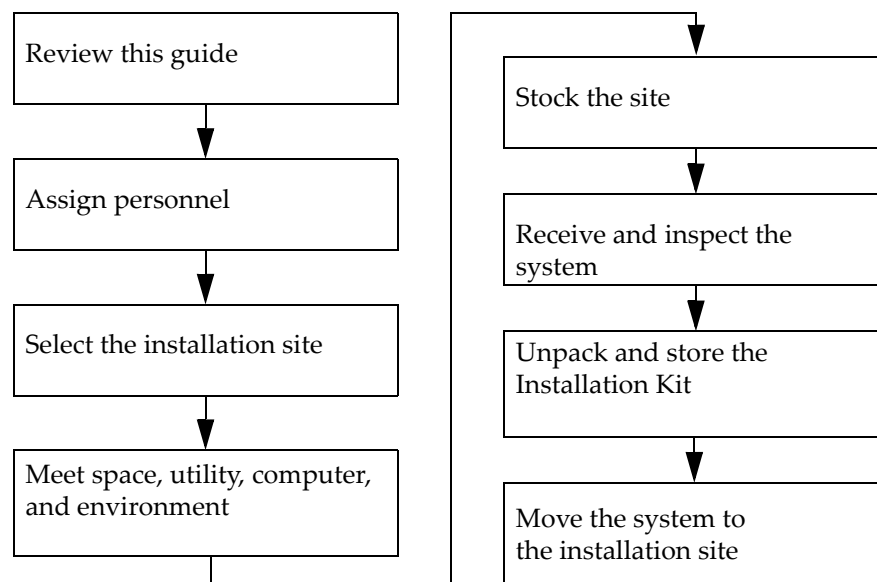
1. Complete the site preparation tasks in this chapter.
2. Fill out the corresponding checklists. See [“Checklists” on page 31](#).
3. Schedule installation before the shipment arrives.
4. Verify with an Applied Biosystems service representative who will contact you by telephone that:
 - All checklists are complete
 - The purchase order is complete
 - You have considered all components and options in preparing the site

Site preparation process

The general site preparation tasks and a suggested sequence for completing the tasks are summarized in [Figure 1, “Site preparation tasks and their suggested sequence”](#), below. The sequence can vary, but always:

- Review this guide first
- Unpack and store the Installation Kit as soon as you receive it

Figure 1 Site preparation tasks and their suggested sequence



Assign personnel

Laboratory safety representative


Applied Biosystems requests that a representative from your laboratory be in the vicinity and available to the Applied Biosystems service representative at all times while the service representative is at your facility. The laboratory safety representative should be familiar with laboratory safety procedures and know the location of all the safety equipment.

Tasks and personnel

[Table 1, Suggested personnel tasks](#), summarizes specific site-preparation tasks and suggests the personnel to accomplish the tasks. Use the table to help schedule and manage the site-preparation process.

Table 1 Suggested personnel tasks

| Personnel | Tasks |
|---|--|
| Site preparation-installation coordinator | <ul style="list-style-type: none"> • Orders required materials. • Chooses the site. • Reviews checklists with applicable personnel, then verifies with the Applied Biosystems service representative that the site is properly prepared. • Schedules the installation and informs personnel of the installation date. • Ensures that the site is clear of unnecessary material on the installation day. • Is available to assist the service representative throughout installation. |
| Laboratory safety representative | <ul style="list-style-type: none"> • Ensures that the required safety practices and equipment are in place. • Is available to assist with unpacking and setup. |
| Laboratory personnel-primary users | <ul style="list-style-type: none"> • Ensures that all customer-provided materials for installation are present at the site. • Primary users (responsible for training other users) are available during the installation, so that they can be trained on the instrument. • Primary users have access to an active email account and networked computer in order to acquire software license during installation. |
| Facilities personnel-primary users | <ul style="list-style-type: none"> • Ensures that installation requirements are met for: <ul style="list-style-type: none"> – Space at the installation site – Building clearances – Temperature and humidity – Waste collection – Electrical supply – Computer – Safety and installation materials • At least two people are available to help the Applied Biosystems service representative move and position the system. |

| Personnel | Tasks |
|---|---|
| Network or IT specialist (if the system will be connected to a network) | <ul style="list-style-type: none"> Ensures that two active, tested local area network (LAN) connections are in place before the scheduled installation date. Ensures that network hardware is compatible with an RJ45-type connector. Is available during installation to connect the system to the network. <p> CAUTION! Do not attempt to connect the system components to the network before the Applied Biosystems service representative arrives.</p> |

Select the site

When deciding where to install the instrument, refer to the following sections for site requirements:

- [“Space requirements” on page 19](#)
- [“Environmental requirements” on page 22](#)
- [“Waste collection requirements” on page 23](#)
- [“Electrical requirements” on page 23](#)
- [“Computer requirements” on page 25](#)
- [“Safety and materials” on page 26](#)

IMPORTANT! The site cannot be designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4). Applied Biosystems does not install, service, or repair Applied Biosystems instruments in areas designated BSL-3 or BSL-4.

Space requirements

System components

The shipment shown in [Figure 2](#) includes the:

- Instrument
- Dell® computer, keyboard and mouse
- Dell® flat screen monitor

Setup requirements

A typical setup is shown in [Figure 2](#) and [Figure 3](#). For details on the space requirements, see “[Required clearances](#)” on page 21.

Figure 2 Setup requirements (not to scale)

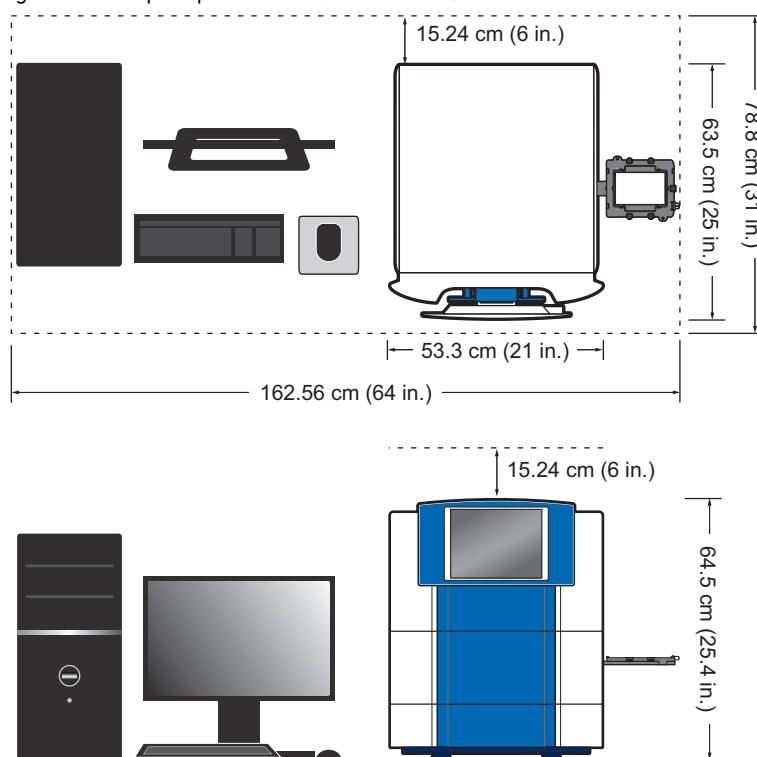
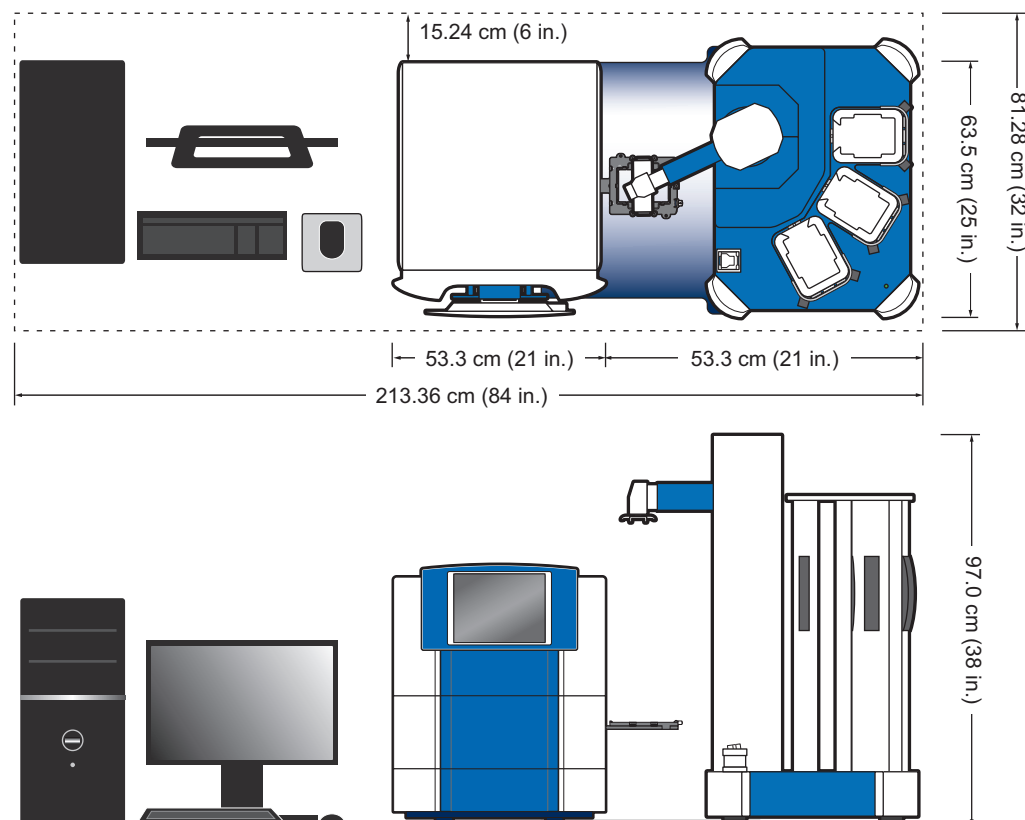


Figure 3 Setup requirements - showing optional robot (not to scale)



Layout requirements

Verify that the installation location:

- Is not adjacent to vibration sources, such as a centrifuge, pump, or compressor. Excessive vibration will affect instrument performance.
- Is not adjacent to electrically noisy devices, such as a refrigeration unit.
- Allows the computer to be within 2 meters (6 feet) of the instrument.
- Allows positioning of the monitor, keyboard, and accessories for proper ergonomics during use.
- Is a stable and level surface. The instrument and robot must be level to function properly.

Dimensions and weights

The table below indicates dimensions and weights of the instrument and associated devices. Ensure that the installation site (floor space and/or bench space) can accommodate the dimensions and support the weights.

Table 2 Dimensions and weights of instrument and associated devices

| Component | Width | Depth | Height | Weight |
|------------------------------|-------------------|--------------------|-------------------|-------------------|
| Instrument | 53.3 cm (21.0 in) | 63.5 cm (25.0 in) | 64.5 cm (25.4 in) | 60.7 kg (134 lbs) |
| Computer (desktop) | 18.7 cm (7.36 in) | 44.5 cm (17.52 in) | 41 cm (16.1 in) | 10.9 kg (24 lbs) |
| Computer (laptop) (optional) | 35.8 cm (14.1 in) | 25.7 cm (10.1 in) | 35.8 cm (14.1 in) | 2.6 kg (5.7 lbs) |
| Computer accessories | | | | |
| • Monitor | 44.7 cm (17.5 in) | 19.3 cm (7.6 in) | 36.6 cm (14.4 in) | 6.9 kg (15.2 lbs) |
| • Keyboard | 44.7 cm (17.5 in) | 15.25 cm (6 in) | 5 cm (2 in) | .09 kg (0.2 lbs) |
| Twister II Robot (optional) | 53.3 cm (21 in) | 77 cm (28 in) | 97 cm (38 in) | 49.9 kg (110 lbs) |

Required clearances

During instrument setup and maintenance, it is necessary to access the back of the instrument. If the back of the instrument faces a wall, it will be necessary to have enough space to rotate the instrument on the bench for access.

Required clearances are:

- Clearance – At least 15.2 cm (6 in) of clearance for ventilation, service access, and cable routing at the back of the instrument. The total width of the required space with the instrument and a desktop computer will be 162.5 cm (64 in) and the total depth will be 81.3 cm (32 in). If a robot is included, the total width will be 213 cm (84 in).
- Vertical clearance – Including instrument height, total vertical space required will be 80 cm (31.4 in). If a robot is included, the total vertical space required will be 112 cm (44 in).

Environmental requirements

Altitude

This instrument is for indoor use only and for altitudes not exceeding 2000 meters (6500 feet) above sea level.

Temperature and humidity requirements

Ensure that the installation site is maintained under the following conditions:

Table 3 Temperature and Humidity requirements

| Condition | Acceptable range |
|-------------|--|
| Temperature | 15 to 30 °C (59 to 86 °F) |
| Humidity | 20 to 80% relative humidity, noncondensing |

Avoid placing the system adjacent to heaters, cooling ducts, or in direct sunlight. Fluctuations between day and night temperatures can cause degradation of performance.

Pollution

The instrument has a Pollution Degree rating of 2. It may be installed in an environment that has nonconductive pollutants only, such as dust particles or wood chips. Typical environments with a Pollution Degree 2 rating are laboratories and sales and commercial areas.

Environmental conditions for storage and transportation

The instrument requires these storage and transport conditions:

- An ambient temperature range of -30°C to +60°C
- A relative humidity range of 20% to 80%

Instrument hot-air exhaust venting

Verify that one of the following conditions exists:

- Facilities personnel have certified that the normal room ventilation system can maintain room temperature if the maximum thermal output of the instrument (2731 Btu/h [800 W]) is vented directly into the room air from the hot-air waste port on the rear panel.
- A suitable venting device such as a fume hood or fume duct is available to vent the hot air exhaust from the instrument space.

Waste collection requirements

Collecting liquid

Dispose of buffer, reagents, and any liquid waste as hazardous waste in compliance with local and national regulations.

For more information, see [“Chemical waste safety” on page 44](#).

Electrical requirements

System electrical requirements

The instrument can accept operating voltages between 100 VAC and 240 VAC with a fluctuation of +/- 10% of the nominal voltage at 50/60 Hz.

Verify that your power supply meets system electrical requirements and perform any required power supply upgrades or changes before the Applied Biosystems service representative arrives to install the instrument.



CAUTION! Do not unpack or plug in any components until the Applied Biosystems service representative has configured the system for the proper operating voltage.

Note: Applied Biosystems recommends that the instrument and computer power receptacle be on an electrical circuit that is not shared with electrically noisy devices or devices that can cause power surges, such as refrigeration units.

[Table 4](#) provides electrical specifications for the instrument and associated devices. For all indicated input voltages, a 15 A circuit is required.

Table 4 Electrical specifications

| Device | Rated voltage (VAC) | Rated Frequency (Hz) | Rated Current (A) | Rated Power (VA) |
|------------------------------|---------------------|----------------------|-------------------|------------------|
| Instrument | 100-240 +/-10% | 50/60 | 12.5 | 950 |
| Computer (desktop) | | | 2.1 | 125 |
| Computer (laptop) (optional) | | | 4.6 | 90 |
| Monitor | | | 1.5 | 65 |
| Twister II Robot (optional) | | | 2.5 | 150 |

Note: The instrument, monitor, desktop computer, robot, and laptop computer will self-adjust for 100v-240v input voltages of 50/60 Hz.

Electrical protective devices

Applied Biosystems recommends several protective devices to protect the system in environments with large voltage and power fluctuations.

Power line regulator Applied Biosystems recommends the use of a 1.5-kVA power line regulator in areas where the supplied power fluctuates in excess of +/- 10% of the normal voltage. Power fluctuations can adversely affect the function of the instrument and computer.

Note: A power line regulator monitors the input current and adjusts the power supplied to the instrument or computer. It does not protect against a power surge or failure.

Uninterruptible power supply (UPS) Applied Biosystems recommends the use of a 1.5-kVA uninterruptible power supply (UPS), especially in areas prone to power failure. Power failures and other events that abruptly terminate the function of the instrument and computer can corrupt data and possibly damage the system.

IMPORTANT! UPSs provide power for a limited time. They are meant to delay the effects of a power outage, not to serve as replacement power sources. In the event of a power loss, power off the instrument and computer unless you expect to regain power within the battery life of the UPS.

Surge protector Applied Biosystems recommends the use of a 10-kVA surge protector (line conditioner) in areas with frequent electrical storms or near devices that are electrically noisy, such as refrigerators, air conditioners, or centrifuges. Short-duration, high-voltage power fluctuations can abruptly terminate the function of, and thereby damage the components of, the computer and the instrument.

Note: A dedicated line and ground between the instrument, computer, and the building's main electrical service can also prevent problems caused by power fluctuations.

Computer requirements

Antivirus software requirements

No antivirus software is provided because customer preferences and network requirements vary. Therefore, you need to install antivirus software of your choice to protect the computer against viruses.



CAUTION! Do not install on the computer additional software other than antivirus software. Changes to the configured software could void the instrument warranty and cause the system to be non-operational.

Network requirements

LAN connection The instrument and computer will be connected to a LAN. Two active, tested LAN connections must be in place before the scheduled installation date. Due to differences in network connections, the Applied Biosystems service representative cannot configure the system to access a specific network.

Network cables The instrument and computer are factory configured for the TCP/IP protocol. The product includes a fast Ethernet adapter (10/100baseT) with an RJ45-type connector and two Ethernet cables that connect the computer and the instrument to the LAN.

IMPORTANT! Do not use a wireless network. A wireless network may interfere with data collection, resulting in data loss.

Printer requirements The instrument and computer can use a network or dedicated printer. The printer and any necessary print drivers must be available before the scheduled installation.

Safety and materials

Safety practices

IMPORTANT! The site must not be designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4). Applied Biosystems does not install, service, or repair Applied Biosystems instruments in areas designated BSL-3 or BSL-4.

IMPORTANT! A safety representative from your facility must ensure that:

- Personnel establish and follow all applicable safety practices and policies to protect laboratory personnel from potential hazards
 - All applicable safety devices and equipment are available at all times
-

Required safety equipment

Your laboratory has specific safety practices and policies designed to protect laboratory personnel from potential hazards that are present. Applied Biosystems expects that you will follow all applicable safety-related procedures at all times.

The following safety protection and equipment must be available at the installation site:

- Protection from any sources of hazardous chemicals, radiation (for example, lasers, radioisotopes, radioactive wastes, and contaminated equipment), and potentially infectious biological material that may be present in the area where the Applied Biosystems service representative will work.
- Appropriate fire extinguisher:
 - You are responsible for providing an appropriate fire extinguisher for use on or near Applied Biosystems equipment.
 - The types and sizes of fire extinguishers shall be suitable for use on electrical and chemical fires as specified in current codes, regulations, and/or standards, and with approval of the Fire Marshall or other authority having jurisdiction.
 - The installation of appropriate fire extinguishers shall be in addition to other fire-protection systems and not as a substitute or alternative to them.
- Eyewash
- Safety shower
- Eye and hand protection
- Adequate ventilation, including vent line/fume hood, if applicable
- Biohazard waste container, if applicable
- First-aid equipment
- Spill cleanup equipment
- Applicable SDSs

Materials for installation

Provide the following materials for the installation:

- Safety glasses, lab coats, chemical-resistant, disposable gloves (powder-free)
- Glassware washing solution
- Lint-free tissues
- Methanol or isopropanol, HPLC-grade or better
- Water, Milli-Q[®] grade
- Three sizes of micropipettors and tips
 - 1 to 10- μ L
 - 10 to 100- μ L
 - 100 to 1,000- μ L
- Mini vortexer, plate adapter, centrifuge, and sample tubes

Materials for routine operation

Additional supplies and consumables are necessary for routine operation of the instrument. Contact the Applied Biosystems sales representative to order these additional supplies.

Use only supplies as specified by Applied Biosystems.

Receiving and inspecting the system

Shipped contents

All shipments include the:

- Instrument
- Sample block
- Heated cover
- Plate adaptor
- Desktop computer (laptop model optional), keyboard, mouse, mouse pad, and monitor
- Calibration reagents
- Instrument verification reagents
- Starter kit
- Software kit
- Power cord
- Array card block (optional)
 - Centrifuge bucket/Clip set (2 types, user selectable)
 - Array card staker
- (Optional) Twister® II Robot with instrument adapter and fixed barcode reader
 - Tall stacks OR short stacks (user selectable)
 - Handheld barcode reader

Shipping list

Verify that the items shown on the shipping list are the same items that you ordered at the time of purchase.

Inspect shipping containers for damage

Carefully inspect the shipping containers and report any damage to the Applied Biosystems service representative. Record any damage or mishandling on the shipping documents.

Unpack and store the Installation Kit

The Installation Kit is boxed separately from the instrument components. When you receive the shipment, unpack the Installation Kit immediately. Store the components as specified in the instructions included with the kit.

IMPORTANT! The installation kit contains reagents that require refrigeration. Upon receiving the shipment, immediately store them as listed in the reagents inserts. Except for the installation kit, do not unpack shipping containers to protect you from liability if any damage occurred during shipping.



CAUTION! Some chemicals used with Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death. Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, work with, or dispose of any chemicals or hazardous materials.

Inspecting shipping containers for damage

Carefully inspect the shipping containers and report any damage to the Applied Biosystems service representative. Record any damage or mishandling on the shipping documents.

Moving the crated instrument to the laboratory

Moving schedule

Before the date of installation:

- Clear the installation site of all unnecessary materials.
- If possible, move the crated instrument from the receiving area to the installation site. Do not uncrate. If possible, move the other shipping containers from the shipping area to the installation site.

Required building clearances

The largest container (crate) included with the shipment contains the instrument. To move the crate to the installation site, verify that the building clearances allow passage of the following crate dimensions:

| Crate dimension | Minimum building clearance |
|-----------------|----------------------------|
| Height | 95 cm (37.4 in) |
| Length | 68.6 cm (27.0 in) |
| Depth | 81.5 cm (32.0 in) |

Instrument weight

The instrument weighs approximately 60.7 kg (134 lbs).

Moving and lifting the instrument



CAUTION! PHYSICAL INJURY HAZARD. The instrument is to be moved and positioned only by the Applied Biosystems service representative. If you decide to lift or move the instrument after it has been installed, do not attempt to lift or move the instrument without the assistance of others, the use of appropriate moving equipment, and proper lifting techniques. Improper lifting can cause painful and permanent back injury. Depending on the weight, moving or lifting an instrument may require two or more persons.



CAUTION! Do not tip the instrument on end. Tipping damages the instrument hardware and electronics.

During installation

After the system is uncrated by your Applied Biosystems Service representative, installation and testing of the instrument and computer takes approximately 4 hours.



CAUTION! While the instrument and computer are being installed, avoid exposure to hazards that may be associated with the installation process.

Operator training

During and/or after installation, the Applied Biosystems service representative reviews data and provides some basic operator training. For additional training and reference information, see the user documents provided with the instrument.

2

Checklists

| | |
|---|----|
| ■ Overview | 32 |
| ■ Personnel checklist | 32 |
| ■ Space and layout checklist | 32 |
| ■ Environmental checklist | 33 |
| ■ Electrical checklist | 33 |
| ■ Computer checklist | 33 |
| ■ Safety checklist | 34 |
| ■ Materials checklist | 34 |
| ■ System receipt and inspection checklist | 35 |
| ■ Moving the crated instrument checklist | 35 |

Overview

Before using the checklists, read all previous sections in this guide.

Use the checklists in this chapter to ensure that you have made all preparations for installing the system. An Applied Biosystems service representative will contact you to verify that all checklists are complete before setting up the installation date.

Personnel checklist

For more information, see [“Assign personnel” on page 16](#).

| Date verified | Designated personnel |
|---------------|---|
| | Site preparation/installation coordinator |
| | Laboratory safety representative |
| | Laboratory personnel: <ul style="list-style-type: none"> To ensure that customer-supplied materials are on hand Primary users to be trained during installation and to subsequently train other users |
| | Facilities personnel: <ul style="list-style-type: none"> To provide environmental, electrical, and computer site-preparation requirements Two people to help the service representative move and position the instrument, if applicable |
| | Network or IT specialist (only if the system will be connected to a network) |

Space and layout checklist

For more information, see [“Space requirements” on page 19](#).

| Date verified | Requirements |
|---------------|---|
| | Location is away from: <ul style="list-style-type: none"> Heating or cooling ducts Direct sunlight Vibration from other instruments or devices Computer workspace allows for proper ergonomics during use. |
| | Location accommodates the dimensions and weights specified in “Dimensions and weights” on page 21 . |
| | Location meets the requirements specified in “Space requirements” on page 19 . |

Environmental checklist

For more information, see [“Environmental requirements” on page 22](#).

| Date verified | Requirement |
|---------------|--|
| | The altitude does not exceed 2000 m (6500 ft) above sea level. |
| | The conditions specified in “Environmental requirements” on page 22 have been met. |
| | Pollution Degree II – Only nonconductive pollutants, if any, are present. |
| | Instrument location is not subject to strong mechanical vibrations. |

Electrical checklist

For more information, see [“Electrical requirements” on page 23](#).

| Date Verified | Requirement |
|---------------|---|
| | The main power supply to the instrument can be immediately disconnected. |
| | Appropriate grounded power receptacles are available (see “Electrical requirements” on page 23). IMPORTANT! Notify your Applied Biosystems Field Service Engineer of any unusual power receptacles or power supply conditions before he/she arrives to do the installation. |

Computer checklist

For more information, see [“Computer requirements” on page 25](#).

| Date verified | Requirement |
|--------------------|---|
| Antivirus Software | |
| | Appropriate antivirus software supplied by the customer is available for loading on the computer. |
| Networking | |
| | Two active, tested LAN connections are available. |
| | Network hardware is compatible with an RJ45-type connector. See “Network requirements” on page 25 . |
| Printer | |
| | If applicable, a network printer or a dedicated printer and necessary print drivers are available. |

Safety checklist

For more information, see [“Safety and materials” on page 26](#).

| Date verified | Requirement |
|---------------|---|
| | The site is not designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4). |
| | Safety practices and policies to protect laboratory personnel from potential hazards are in place and are followed. |
| | Protection from any sources of hazardous chemicals, radiation (for example, lasers, radioisotopes, radioactive wastes, and contaminated equipment), and potentially infectious biological material is in place. |
| | Appropriate fire extinguisher |
| | Eye and hand protection |
| | Eyewash |
| | Safety shower |
| | Vent lines/fume hood, if applicable |
| | Biohazard waste container, if applicable |
| | First-aid equipment |
| | Spill cleanup equipment |
| | SDSs readily available |

Materials checklist

For more information, see [“Safety and materials” on page 26](#).

| Date Verified | Requirement |
|------------------------------------|---|
| Materials for General Installation | |
| | Safety glasses and lab coats |
| | Chemical-resistant disposable gloves (powder free) |
| | Glassware washing solution |
| | Lint-free tissues |
| | Methanol or isopropanol, HPLC-grade or better |
| | Water, Milli-Q® grade |
| | Three sizes of micropipettors and tips: <ul style="list-style-type: none"> • 1- to 10-µL • 10- to 100-µL • 100- to 1000-µL |
| | A mini vortexer, plate adapter, centrifuge, and sample tubes |

| Date Verified | Requirement |
|---------------------------------|--|
| Materials for Routine Operation | |
| | Materials for routine operation after the installation are available or have been ordered (see “Materials for routine operation” on page 27). |


System receipt and inspection checklist

For more information, see [“Receiving and inspecting the system” on page 28](#).

| Date verified | Action |
|---------------|--|
| | Verified that items on the packing list are those that were ordered. Otherwise, reported to the Applied Biosystems service representative discrepancies in the packing list. |
| | Opened and stored the Installation Kit components as specified in the kit operating instructions. |
| | Received the system and inspected the shipping containers for mishandling or damage. IMPORTANT! Do not open any shipping containers except for the Installation Kit. |
| | Reported to the Applied Biosystems service representative any signs of mishandling or damage to the shipping container. |

Moving the crated instrument checklist

For more information, see [“Moving the crated instrument to the laboratory” on page 29](#).

| Date Verified | Item |
|---------------|---|
| | The measured building clearances can accommodate the instrument crate dimensions (see “Required building clearances” on page 29). If the crate dimensions exceed building clearances, contact the Applied Biosystems service representative. Do not unpack the crate without authorization. |
| | If possible, move all the crated equipment, excluding the crated instrument, to the laboratory before the date of the scheduled installation.  WARNING! PHYSICAL INJURY HAZARD. Do not attempt to lift or move any boxed or crated items unless you have received related training. Incorrect lifting can cause painful and sometimes permanent back injury. Use proper lifting techniques when lifting or moving items. |
| | Cleared the installation site of all unnecessary materials. |



Safety and EMC Compliance Information







| | |
|--|----|
| ■ Symbols on instruments | 38 |
| ■ General instrument safety | 41 |
| ■ Chemical safety | 42 |
| ■ Chemical waste safety | 44 |
| ■ Electrical safety | 45 |
| ■ Physical hazard safety | 46 |
| ■ Biological hazard safety | 47 |
| ■ Workstation safety | 47 |
| ■ Safety and electromagnetic compatibility (EMC) standards | 48 |



Symbols on instruments




Electrical symbols on instruments





The following table describes the electrical symbols that may be displayed on Applied Biosystems instruments.

| Symbol | Description |
|--|--|
|  | Indicates the On position of the main power switch. |
|  | Indicates the Off position of the main power switch. |
|  | Indicates a terminal that may be connected to the signal ground reference of another instrument. This is not a protected ground terminal. |
|  | Indicates a protective grounding terminal that must be connected to earth ground before any other electrical connections are made to the instrument. |
|  | Indicates a terminal that can receive or supply alternating current or voltage. |
|  | Indicates that the device receives or supplies direct current or voltage. |

Safety symbols


The following table describes the safety symbols that may be displayed on Applied Biosystems instruments. Each symbol may appear by itself or with text that explains the relevant hazard. These safety symbols may also appear next to DANGERS, WARNINGS, and CAUTIONS that occur in the text of this and other product-support documents.

| Symbol | Description | Français |
|---|--|---|
|  | Caution, risk of danger. Consult the user documentation for additional information and instructions, and proceed with appropriate caution. | ATTENTION! Lire les fiches techniques de sûreté et d'utilisation de matériels avant toute manipulation de produits. Procéder délicatement. |
|  | Caution, risk of electric shock. | ATTENTION! Risque de choc électrique. |
|  | Caution, hot surface. | ATTENTION! Surface brûlante. |

| Symbol | Description | Français |
|--|-----------------------------|--|
|   | Caution, moving parts. | ATTENTION! Pièces en mouvement, risque de pincement et/ou d'écrasement. |
|  | Caution, laser radiation. | ATTENTION! Radiation Laser. |
|  | Caution, ultraviolet light. | ATTENTION! Emission ultra violet. |

Environmental symbols on instruments

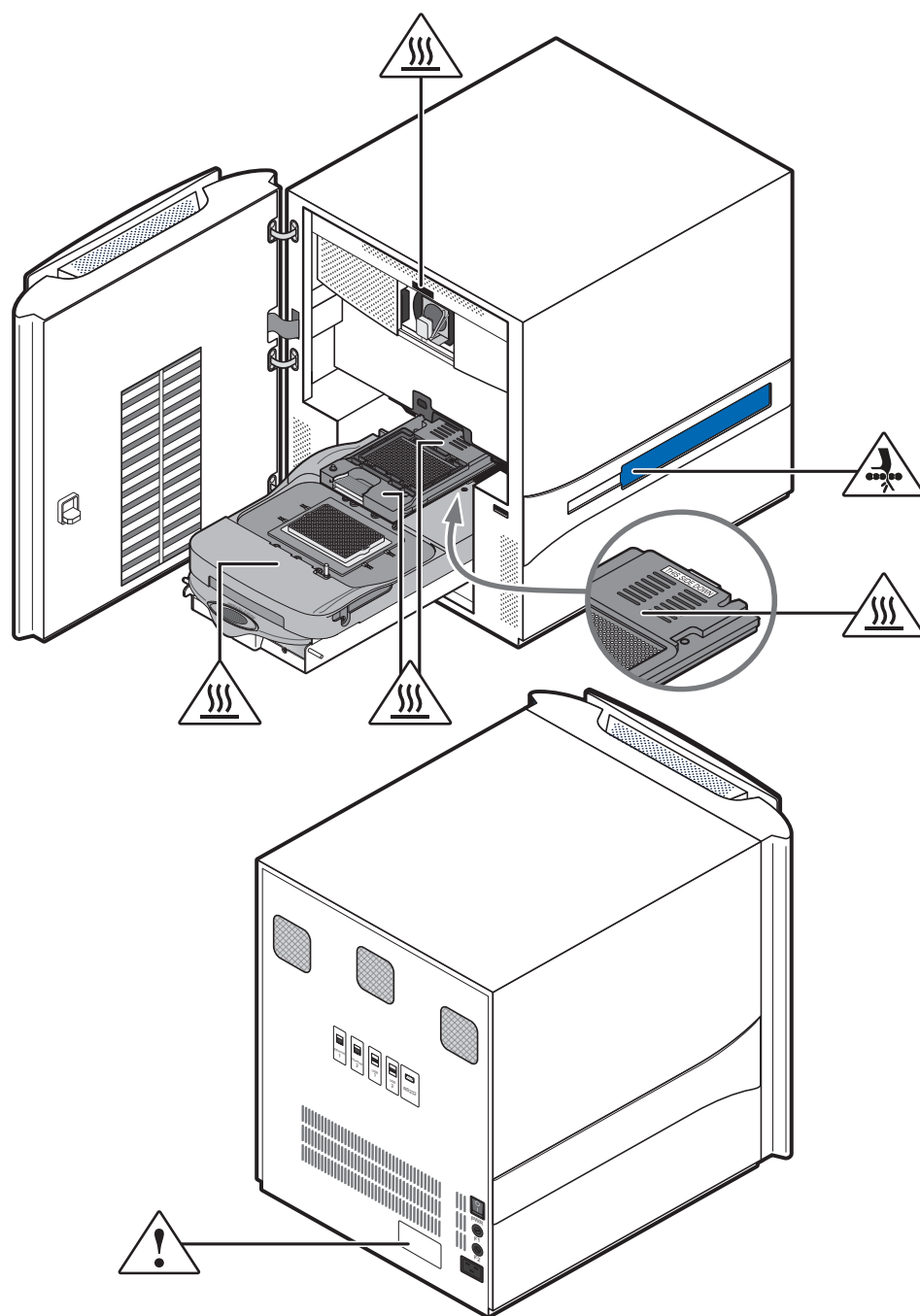
The following symbol applies to all Applied Biosystems electrical and electronic products placed on the European market after August 13, 2005.

| Symbol | Description |
|---|---|
|  | <p>Do not dispose of this product as unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provisions to reduce the environmental impact of waste electrical and electronic equipment (WEEE).</p> <p>European Union customers: Call your local Applied Biosystems Customer Service office for equipment pick-up and recycling. See www.appliedbiosystems.com for a list of customer service offices in the European Union.</p> |



Location of safety symbols

The instrument contains safety symbol warnings at the locations shown below:



General instrument safety



WARNING! Physical Injury Hazard. Using the instrument in a manner not specified by Applied Biosystems may result in personal injury or damage to the instrument.

Moving and lifting the instrument



CAUTION! Physical Injury Hazard. The instrument is to be moved and positioned only by the personnel or vendor specified in the applicable site preparation guide. If you decide to lift or move the instrument after it has been installed, do not attempt to lift or move the instrument without the assistance of others, the use of appropriate moving equipment, and proper lifting techniques. Improper lifting can cause painful and permanent back injury. Depending on the weight, moving or lifting an instrument may require two or more persons.

Moving and lifting stand-alone computers and monitors



CAUTION! Do not attempt to lift or move the computer or the monitor without the assistance of others. Depending on the weight of the computer and/or the monitor, moving them may require two or more people.

Things to consider before lifting the computer and/or the monitor:

- Make sure that you have a secure, comfortable grip on the computer or the monitor when lifting.
- Make sure that the path from where the object is to where it is being moved is clear of obstructions.
- Do not lift an object and twist your torso at the same time.
- Keep your spine in a good neutral position while lifting with your legs.
- Participants should coordinate lift and move intentions with each other before lifting and carrying.
- Instead of lifting the object from the packing box, carefully tilt the box on its side and hold it stationary while someone slides the contents out of the box.

Operating the instrument

Ensure that anyone who operates the instrument has:

- Received instructions in both general safety practices for laboratories and specific safety practices for the instrument.
- Read and understood all applicable Safety Data Sheets (SDSs). See [“Obtain an SDS” on page 42.](#)



Cleaning or decontaminating the instrument



CAUTION! Before using a cleaning or decontamination method other than those recommended by the manufacturer, verify with the manufacturer that the proposed method will not damage the equipment.

Chemical safety

Chemical hazard warning



WARNING! CHEMICAL HAZARD. Before handling any chemicals, refer to the Safety Data Sheet (SDS) provided by the manufacturer, and observe all relevant precautions.



WARNING! CHEMICAL HAZARD. All chemicals in the instrument, including liquid in the lines, are potentially hazardous. Always determine what chemicals have been used in the instrument before changing reagents or instrument components. Wear appropriate eyewear, protective clothing, and gloves when working on the instrument.

About SDSs

Chemical manufacturers supply current Safety Data Sheets (SDSs) with shipments of hazardous chemicals to new customers. They also provide SDSs with the first shipment of a hazardous chemical to a customer after an SDS has been updated. SDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new SDS packaged with a hazardous chemical, be sure to replace the appropriate SDS in your files.

Obtain an SDS

The Safety Data Sheet (SDS) for any chemical supplied by Applied Biosystems is available to you free 24 hours a day. To obtain an SDS:

1. Go to www.appliedbiosystems.com, click **Support**, then select **SDS**.
2. In the Keyword Search field of the SDS Search page:
 - a. Type in the chemical name, part number, or other information that you expect to appear in the SDS of interest.
 - b. Select the language of your choice.
 - c. Click **Search**.

A list of relevant SDS documents appears.

3. Find the document of interest, right-click the document title, then select any of the following:

- **Open** – To view the document
- **Save Target As** – To download a PDF version of the document to a destination that you choose
- **Print Target** – To print the document

Note: For the SDSs of chemicals not distributed by Applied Biosystems, contact the chemical manufacturer.

4. To have a copy of an SDS sent by email, send your request to:
SDS_Inquiry_CCRM@appliedbiosystems.com
5. To have a copy of an SDS sent by fax to the US or Canada, send your fax request to 650-442-2252.
6. To have a copy of an SDS sent by fax to a region outside of the US or Canada, scroll to the bottom of your SDS Search results web page and click **regional office**.

From the Country/Region drop-down menu, select your location.

The contact phone number and the fax request number for your region appear.

Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See “Obtain an SDS” on page 42.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the SDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the SDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer’s cleanup procedures as recommended in the SDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.



Chemical waste safety

Chemical waste hazard



CAUTION! HAZARDOUS WASTE. Refer to Safety Data Sheets and local regulations for handling and disposal.



WARNING! CHEMICAL WASTE HAZARD. Wastes produced by Applied Biosystems instruments are potentially hazardous and can cause injury, illness, or death.



WARNING! CHEMICAL STORAGE HAZARD. Never collect or store waste in a glass container because of the risk of breaking or shattering. Reagent and waste bottles can crack and leak. Each waste bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position. Wear appropriate eyewear, clothing, and gloves when handling reagent and waste bottles.

Chemical waste safety guidelines

To minimize the hazards of chemical waste:

- Read and understand the Safety Data Sheets (SDSs) provided by the manufacturers of the chemicals in the waste container before you store, handle, or dispose of chemical waste.
- Provide primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing). For additional safety guidelines, consult the SDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the SDS.
- Handle chemical wastes in a fume hood.
- After emptying a waste container, seal it with the cap provided.
- Dispose of the contents of the waste tray and waste bottle in accordance with good laboratory practices and local, state/provincial, or national environmental and health regulations.

Waste disposal

If potentially hazardous waste is generated when you operate the instrument, you must:

- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure the health and safety of all personnel in your laboratory.
- Ensure that the instrument waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.

IMPORTANT! Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.

Electrical safety



DANGER! ELECTRICAL SHOCK HAZARD. Severe electrical shock can result from operating the instrument without its instrument panels in place. Do not remove instrument panels. High-voltage contacts are exposed when instrument panels are removed from the instrument.

Power



DANGER! ELECTRICAL HAZARD. Grounding circuit continuity is vital for the safe operation of equipment. Never operate equipment with the grounding conductor disconnected.



DANGER! ELECTRICAL HAZARD. Use properly configured and approved line cords for the voltage supply in your facility.



DANGER! ELECTRICAL HAZARD. Plug the system into a properly grounded receptacle with adequate current capacity.

Overvoltage rating

The instrument and associated devices have an installation (overvoltage) category of II, and are classified as portable equipment.



Fuses



WARNING! FIRE HAZARD. Improper fuses or high-voltage supply can damage the instrument wiring system and cause a fire. Before turning on the instrument, verify that the fuses are properly installed and that the instrument voltage matches the power supply in your laboratory.



WARNING! FIRE HAZARD. For continued protection against the risk of fire, replace fuses only with fuses of the type and rating specified for the instrument.

Bar code scanner laser safety

Laser
classification

The bar code scanners included with the instrument are categorized as Class 2 (II) lasers.

Laser safety
requirements

Class 2 (II) lasers are low-power, visible-light lasers that can damage the eyes. Never look directly into the laser beam. The scanner is designed to prevent human access to harmful levels of laser light during normal operation, user maintenance, or during prescribed service operations.



WARNING! LASER HAZARD. Class 2 (II) lasers can cause damage to eyes. Avoid looking into a Class 2 (II) laser beam or pointing a Class 2 (II) laser beam into another person's eyes.

Physical hazard safety

Ultraviolet light



WARNING! ULTRAVIOLET LIGHT HAZARD. Looking directly at a UV light source can cause serious eye damage. Never look directly at a UV light source and always prevent others from UV exposure. Follow the manufacturer's recommendations for appropriate protective eyewear and clothing.

Moving parts



WARNING! PHYSICAL INJURY HAZARD. Moving parts can crush and cut. Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing the instrument.

Biological hazard safety

General biohazard



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective equipment, which includes but is not limited to: protective eyewear, face shield, clothing/lab coat, and gloves. All work should be conducted in properly equipped facilities using the appropriate safety equipment (for example, physical containment devices). Individuals should be trained according to applicable regulatory and company/institution requirements before working with potentially infectious materials. Read and follow the applicable guidelines and/or regulatory requirements in the following:

- U.S. Department of Health and Human Services guidelines published in *Biosafety in Microbiological and Biomedical Laboratories* (stock no. 017-040-00547-4;)
<http://www.cdc.gov/biosafety/publications/index.htm>
- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html).
- Your company's/institution's Biosafety Program protocols for working with/handling potentially infectious materials.

Additional information about biohazard guidelines is available at:

www.cdc.gov

Workstation safety

Correct ergonomic configuration of your workstation can reduce or prevent effects such as fatigue, pain, and strain. Minimize or eliminate these effects by configuring your workstation to promote neutral or relaxed working positions.



WARNING! MUSCULOSKELETAL AND REPETITIVE MOTION HAZARD. These hazards are caused by potential risk factors that include but are not limited to repetitive motion, awkward posture, forceful exertion, holding static unhealthy positions, contact pressure, and other workstation environmental factors.

To minimize musculoskeletal and repetitive motion risks:

- Use equipment that comfortably supports you in neutral working positions and allows adequate accessibility to the keyboard, monitor, and mouse.
- Position the keyboard, mouse, and monitor to promote relaxed body and head postures.



Safety and electromagnetic compatibility (EMC) standards

This section provides information on:

- [U.S. and Canadian safety standards](#)
- [Canadian EMC standard](#)
- [European safety and EMC standards](#)
- [Australia and New Zealand EMC standards](#)

U.S. and Canadian safety standards



The instrument has been tested to and complies with standard:

UL 61010-1:2nd Edition/CSA C22.2 No. 61010-1, "Safety Requirements for Electrical Equipment for Measurement, Control, and Laboratory Use, Part 1: General Requirements."

UL 61010-2-010, "Particular Requirements for Laboratory Equipment for the Heating of Materials."

Canadian EMC standard

This instrument has been tested to and complies with standard:

ICES-001, Issue 3: "Industrial, Scientific, and Medical Radio Frequency Generators." Cet appareil numérique de la classe B est conforme à la norme NMB-001 du Canada.

European safety and EMC standards

Safety

This instrument meets European requirements for safety (Low Voltage Directive 2006/95/EC). This instrument has been tested to and complies with standards:

EN 61010-1:2001, "Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use, Part 1: General Requirements."

EN 61010-2-010:2003, "Particular Requirements for Laboratory Equipment for the Heating of Materials."

EN 61010-2-081:2002+A1:2003, "Particular Requirements for Automatic and Semi-Automatic Laboratory Equipment for Analysis and Other Purposes."

EMC

The ViiA™ 7 Real-Time PCR System meets European requirements for emission and immunity (EMC Directive 2004/108/EC).

EN 61326-1:2006 "Electrical equipment for measurement, control and laboratory use-Part 1 General EMC requirements." (Group 1, Class B)

Australia and New Zealand EMC standards



This instrument has been tested to and complies with standard AS/NZS 2064, "Limits and Methods Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radio-frequency Equipment."





Appendix A
Safety and electromagnetic compatibility (EMC) standards

A

antivirus software 25

B

biohazardous waste, handling 45

biosafety levels 18

building clearances, required 29

C

CAUTION, description 8

checklists

computer 33

electrical 33

environmental 33

inspection 35

layout 32

materials 34

moving the instrument 35

personnel assignments 32

safety 34

site 32

site preparation 31

space 32

system receipt 35

chemical safety 42, 43

chemical waste safety 44

clearances

building 29

required 21

system components 21

computer checklists 33

consumables, ordering 27

conventions

IMPORTANT! 7

Notes 7

safety 8

user attention words 7

crate

See also moving the crated system

dimensions 29

reporting shipping damage or tipping 29

customer training, information 30

D

damage, reporting 29

DANGER, description 8

E

electrical checklist 33

electrical protective devices 24

electrical safety 45

electrical specifications 23

electromagnetic compatibility standards. *See* EMC standards

EMC standards 48

environmental checklist 33

environmental requirements 22

ergonomics, safety 47

F

facilities personnel, tasks 16

field service, installation tasks 30

fire extinguisher 26

G

General 47

guidelines

chemical safety 43

chemical waste disposal 44

chemical waste safety 44

H

hazard icons. *See* safety symbols, on instruments

hazard symbols. *See* safety symbols, on instruments

hazards. *See* safety

I

IMPORTANT, description 8
 inspection checklist 35
 installation
 conditions for postponing 14
 field service tasks 30
 overview 30
 scheduling 14
 time required 30
 installation category 45
 installation coordinator
 personnel assignments 32, 34
 instrument
 basic operator training 30
 components shipped 19, 28
 instrument operation, safety 41
 IT specialist, tasks 17

L

laboratory layout, typical 20
 laboratory personnel, tasks 16
 laboratory safety representative 16
 laser classification 46
 laser safety
 bar code scanner 46
 requirements 46
 layout checklist 32
 lifting, safety 29
 line conditioner, requirements 24
 installation site. *See* laboratory layout

M

materials
 checklist 34
 for routine operation 27
 moving and lifting, safety 41
 moving parts, safety 46
 moving the crated system
 building clearances 29
 no lifting caution 29
 safety 29
 weight 29
 moving the instrument checklist 35

N

network (IT) specialist, tasks 17

network specialist, tasks 17

O

overvoltage category (rating) 45

P

personnel assignments 16, 32
 personnel checklist 32
 physical hazard safety 46
 power line regulator 24
 protective devices, electrical 24

R

radioactive waste, handling 45
 regulator, power line 24
 repetitive motion, safety 47
 reporting, shipping damage 29
 requirements
 antivirus software 25
 component clearances and positioning 21
 environmental 22
 installation 30
 no tipping 29
 physical clearances 21
 quick disconnect 18
 space 19, 21
 temperature and humidity 22

S

safety
 bar code scanner 46
 before operating the instrument 41
 biological hazards 47
 checklist 34
 chemical 42
 chemical waste 44
 conventions 8
 electrical 45
 ergonomic 47
 guidelines 43, 44, 45
 instrument operation 41
 laser 46
 moving and lifting 29
 moving and lifting instrument 41
 moving parts 46
 moving/lifting 41

- physical hazard 46
- repetitive motion 47
- standards 48
- ultraviolet light 46
- workstation 47
- safety devices, availability 26
- safety equipment, availability 26
- safety information
 - electrical quick disconnect 18
 - lifting 29
- safety labels, on instruments 10
- safety practices, required 26
- safety standards 48
- safety symbols, on instruments 38
- SDSs
 - description 42
 - obtaining 11
- shipping crate. *See* Crate
- shipping damage, reporting 29
- shipping list, verifying 19, 28
- shipping, reporting damage 29
- site
 - checklist 32
- site preparation
 - See also* requirements
 - checklists 31
 - flowchart 15
 - overview 14
 - schedule 14
 - tasks 16
 - tasks overview 14
- site preparation process, overview 15
- space checklist 32
- space requirements 19, 21
- specifications
 - electrical 23
- standards
 - EMC 48
 - safety 48
- surge protector, requirements 24
- system
 - building clearances 29
 - components 19
 - dimensions 21
 - requirements 18
 - weight 21
- system receipt checklist 35

T

- tasks
 - facilities personnel 16
 - IT specialist 17
 - laboratory personnel 16
 - laboratory safety representative 16
 - network specialist 17
- temperature and humidity, acceptable range 22
- Test Standards Kit, unpacking and storing 28
- training
 - information on 11
 - operator 30

U

- ultraviolet light, safety 46
- uninterruptible power supply, requirements 24
- unpacking, test standards kit 28
- UPS, requirements 24
- user attention words, described 7

W

- WARNINGS
 - chemical hazards 28
 - lifting 29
 - tipping the instrument 29
- waste disposal, guidelines 45
- waste profiles, description 45
- workstation safety 47

Part Number 4445302 Rev B



Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA
Phone 760.603.7200
www.lifetechnologies.com

Technical Resources and Support

For the latest technical resources and support information
for all locations, please refer to our Web site at
www.appliedbiosystems.com/support