# **applied**biosystems

# StepOne<sup>™</sup> and StepOnePlus<sup>™</sup> Real-Time PCR Systems SITE PREPARATION GUIDE

Publication Number 4376768

Revision E



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**

Table 1 Revision history of Pub. no. 4376768

Revision	Date	Description
Е		Minor update in environmental requirements. Update in logos and covers.

#### Limited Use Label License No. 474: Real-Time PCR System for Research Use Only

Notice to Purchaser: The purchase of this instrument conveys to the purchaser the limited, non-transferable right to use the purchased instrument only, under intellectual property rights that are owned and/or controlled by Life Technologies and relate specifically to the instrument. Purchase of the instrument includes the right to use the instrument for internal research and to perform services (including the right to report the results of services for a fee) by the purchaser only, but does not convey rights to use any other products, reagents, assays or methods such as the 5' nuclease assay process. The sale of this instrument is expressly conditioned on the purchaser not reselling, repackaging, or distributing this instrument, or any of its components, and no such rights are conveyed expressly, by implication, or by estoppel. For information on obtaining additional rights, please contact **outlicensing@lifetech.com** or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

#### **Corporate entity**

Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

#### **TRADEMARKS**

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2016 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

		<b>F</b> ~	
$\boldsymbol{P}$	$r \omega$	12	റമ
	rei	ш	-

	How to Use This Guide	5
	How to Obtain More Information	6
	How to Obtain Support	7
	Safety Conventions Used in This Document	8
	Symbols on Instruments	9
	Safety Labels on Instruments	10
	General Instrument Safety	11
	Chemical Safety	12
	Chemical Waste Safety	14
	Electrical Safety	15
	Biological Hazard Safety	15
	LED Safety Requirements	16
	Workstation Safety	16
	Safety and Electromagnetic Compatibility (EMC) Standards	17
Chantar 1	Cita Dranguation Tools	
Chapter 1	Site Preparation Tasks	
Chapter 1	Site Preparation Tasks  Overview	. 1-2
Chapter 1	·	
Chapter 1	Overview	. 1-3
Chapter 1	Overview	. 1-3 . 1-4
Chapter 1	Overview Assigning Personnel Selecting the Site	. 1-3 . 1-4 . 1-4
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements	. 1-3 . 1-4 . 1-4 . 1-7
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements	. 1-3 . 1-4 . 1-4 . 1-7
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements Electrical Requirements	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements Electrical Requirements Computer Requirements	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8 . 1-9
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements Electrical Requirements Computer Requirements Network Requirements	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8 . 1-9 1-10
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements Electrical Requirements Computer Requirements Network Requirements Safety and Materials	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8 . 1-9 . 1-10
Chapter 1	Overview  Assigning Personnel Selecting the Site Space Requirements Environmental Requirements Ventilation Requirements Electrical Requirements Computer Requirements Network Requirements Safety and Materials Receiving and Inspecting the System	. 1-3 . 1-4 . 1-7 . 1-8 . 1-8 . 1-9 . 1-10 1-11

# Chapter 2 Checklists

Overview	2-2
Personnel Checklist	2-2
Space and Layout Checklist	2-3
Environmental Checklist	2-3
Ventilation Checklist	2-3
Electrical Checklist	2-4
Computer and Networking Checklist	2-4
Safety Checklist	2-5
Materials Checklist	2-6
System Receipt and Inspection Checklist	2-7
Moving the Crated Instrument Checklist	2-7

# Index

# **Preface**

### How to Use This Guide

# Purpose of This Guide

This guide helps you prepare to receive and install the Applied Biosystems StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System.

#### Audience

This guide is intended for personnel who will schedule, manage, and perform the tasks required to prepare your site for installation of the StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems.

#### **Assumptions**

This guide assumes that you have:

- Familiarity with Microsoft Windows® XP operating system.
- Familiarity with the Internet and internet browsers.
- Knowledge of techniques for handling DNA samples and preparing them for PCR.
- An understanding of data storage, file transfer, and copying and pasting.
- Networking experience, if you plan to integrate the StepOne<sup>TM</sup> system into your existing laboratory data flow.

#### **Text Conventions**

This guide uses the following conventions:

- **Bold** text indicates user action. For example:
  - Type **0**, then press **Enter** for each of the remaining fields.
- *Italic* text indicates new or important words and is also used for emphasis. For example:
  - Before analyzing, always prepare fresh matrix.
- A right arrow symbol (▶) separates successive commands you select from a drop-down or shortcut menu. For example:
  - Select File ▶ Open ▶ Spot Set.

# User Attention Words

Two user attention words appear in the 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.

### How to Obtain More Information

### Related Documentation

The following documentation ships with the StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems:

Document	PN
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Getting Started Guide for Genotyping Experiments	4376786
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Getting Started Guide for Presence/Absence Experiments	4376787
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> ( $\Delta\Delta C_T$ ) Experiments	4376785
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Getting Started Guide for Standard Curve Experiments	4376784
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Installation, Maintenance, and Networking Guide	4376782
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Installation Quick Reference Card	4376783
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Site Preparation Guide	4376768
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR Systems Software Help	_

The following supporting StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems documentation is available for purchase:

Document	PN
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR System Installation Performance Verification Protocol	4376791
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR System Installation Qualification-Operation Qualification Protocol	4376790
Applied Biosystems StepOne <sup>™</sup> and StepOnePlus <sup>™</sup> Real-Time PCR System Planned Maintenance Protocol	4376788

**Note:** For additional documentation, see "How to Obtain Support" on page 7.

### Obtaining Information from the Help System

The StepOne<sup>™</sup> Software has a Help system that describes how to use each feature of the user interface. Access the Help system by doing one of the following:

- If displayed, click ② in the toolbar of the StepOne<sup>™</sup> Software window
- Select Help ▶ Contents and Index
- Press F1

You can use the Help system to find topics of interest by:

- Reviewing the table of contents
- Searching for a specific topic
- Searching an alphabetized index

#### Send Us Your Comments

We welcome 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 **http://www.appliedbiosystems.com**, then click the link for **Support**. (See "How to Obtain Support" below).

# How to Obtain Support

For the latest services and support information for all locations, go to <a href="http://www.appliedbiosystems.com">http://www.appliedbiosystems.com</a>, 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 user documents, MSDSs, certificates of analysis, and other related documents
- · 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 Technical Support and Sales facilities.

# Safety Conventions Used in This Document

### Safety Alert Words

Four safety alert words appear in the 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 limited to the most extreme situations.

Except for IMPORTANTs, each safety alert word in the 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 the instruments (see "Safety Symbols" on page 9).

#### **Examples**

**IMPORTANT!** You must create a separate sample entry spreadsheet for each 96-well plate.

**CAUTION** CHEMICAL HAZARD. TaqMan Universal PCR Master Mix may cause eye and skin irritation. Exposure may cause discomfort if swallowed or inhaled. Read the MSDS, and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

WARNING PHYSICAL INJURY HAZARD. During instrument operation, the heated cover and sample block can reach temperatures in excess of 100 °C.

**DANGER** ELECTRICAL HAZARD. Grounding circuit continuity is vital for the safe operation. Never operate the instrument with the grounding conductor disconnected.

# Symbols on Instruments

### Electrical Symbols on Instruments

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

Symbol	Description	Symbol	
	Indicates the <b>On</b> position of the main power switch.	=	Indic conn refere This
	Indicates the <b>Off</b> position of the main power switch.		termi
			Indic termi
0	Indicates a standby switch by which the instrument is switched on to the <b>Standby</b> condition. Hazardous voltage may be	~	to ea elect to the
	present if this switch is on standby.		Indic recei curre
Ф	Indicates the <b>On/Off</b> position of a push-push main power switch.	=	Indic recei direc

### Description cates a terminal that may be nected to the signal ground rence of another instrument. is not a protected ground cates a protective grounding ninal that must be connected arth ground before any other trical connections are made e instrument. cates a terminal that can eive or supply alternating ent or voltage. cates a terminal that can eive or supply alternating or ct current or voltage.

### Safety Symbols

The following table describes the safety symbols that may be displayed on the instruments. Each symbol may appear by itself or with text that explains the relevant hazard (see "Safety Labels on Instruments" on page 10). 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	S
<u></u>	Indicates that you should consult the manual for further information and to proceed with appropriate caution.	
<u> </u>	Indicates the presence of a hot surface or other high-temperature hazard and to proceed with appropriate caution.	2

Symbol	Description
	Indicates the presence of moving parts and to proceed with appropriate caution.
<u></u>	Indicates the presence of an electrical shock hazard and to proceed with appropriate caution.
*	Indicates the presence of a laser inside the instrument and to proceed with appropriate caution.

### Environmental Symbols on Instruments

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

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

# Safety Labels on Instruments

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

English	Français
<b>CAUTION</b> Hazardous chemicals. Read the Material Safety Data Sheets (MSDSs) before handling.	ATTENTION Produits chimiques dangeureux. Lire les fiches techniques de sûreté de matériels avant la manipulation des produits.
CAUTION Hazardous waste. Refer to MSDS(s) and local regulations for handling and disposal.	ATTENTION Déchets dangereux. Lire les fiches techniques de sûreté de matériels et la régulation locale associées à la manipulation et l'élimination des déchets.
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 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 de Thermo Fisher Scientific.
CAUTION Moving parts.	ATTENTION Parties mobiles.
<b>DANGER</b> Class 3B (III) visible and/or invisible LED radiation present when open and interlocks defeated. Avoid exposure to beam.	DANGER Rayonnement visible ou invisible d'un faisceau LED de Classe 3B, (III) en cas d'ouverture et de neutralisation des dispositifs de sécurité. Evitez toute exposition au faisceau.

# Location of Warning

The StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> instruments contain a warning label at the location shown below:



# **General Instrument Safety**

WARNING PHYSICAL INJURY HAZARD. Using the instrument in a manner not specified by Thermo Fisher Scientific may result in personal injury or damage to the instrument.

Moving and Lifting the Instrument 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 Standalone Computers and Monitors **WARNING** 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 Material Safety Data Sheets (MSDSs). See "About MSDSs" on page 13.

### 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 Material Safety Data Sheet (MSDS) provided by the manufacturer, and observe all relevant precautions.

**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 Safety Guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. (See "About MSDSs" on page 13.)
- 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 MSDS.
- 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 MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

#### **About MSDSs**

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

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

# Obtaining MSDSs

The MSDS for any chemical supplied by us is available to you free 24 hours a day. To obtain MSDSs:

- 1. Go to https://docs.appliedbiosystems.com/msdssearch.html
- 2. In the Search field of the MSDS Search page:
  - a. Type in the chemical name, part number, or other information that you expect to appear in the MSDS of interest.
  - b. Select the language of your choice.
  - c. Click Search.
- 3. To view, download, or print the document of interest:
  - a. Right-click the document title.
  - b. Select:
    - 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
- 4. To have a copy of an MSDS sent by fax or e-mail, in the Search Results page:
  - a. Select Fax or Email below the document title.
  - b. Click **RETRIEVE DOCUMENTS** at the end of the document list.
  - c. Enter the required information.
  - d Click View/Deliver Selected Documents Now

**Note:** For the MSDSs of chemicals not distributed by us, contact the chemical manufacturer.

# **Chemical Waste Safety**

#### Chemical Waste Hazard

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

WARNING CHEMICAL WASTE HAZARD. Wastes produced by the 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 Material Safety Data Sheets (MSDSs) 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 MSDS.
- 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 MSDS.
- 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 StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems without their instrument panels in place. Do not remove instrument panels. High-voltage contacts are exposed when instrument panels are removed from the instrument.

#### **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.

#### **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 instrument into a properly grounded receptacle with adequate current capacity.

### Overvoltage Rating

The StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems has an installation (overvoltage) category of II, and is classified as portable equipment

# **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://bmbl.od.nih.gov)

- Occupational Safety and Health Standards, Bloodborne Pathogens (29 CFR§1910.1030; http://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:

http://www.cdc.gov

# **LED Safety Requirements**

To ensure safe LED operation:

- The system must be maintained by a Thermo Fisher Scientific Technical Representative.
- All instrument panels must be in place on the instrument while the instrument is operating. When all panels are installed, there is no detectable radiation present. If any panel is removed when the LED is operating (during service with safety interlocks disabled), you may be exposed to LED emissions in excess of the Class **3B** rating.
- Do not remove safety labels or disable safety interlocks.

# **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 vour workstation to promote neutral or relaxed working positions.

### CAUTION 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

U.S. and Canadian Safety Standards The StepOne<sup>TM</sup> and StepOnePlus<sup>TM</sup> systems has been tested to and complies with standard:



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

FDA "Radition Control for Health and Safety Act of 1968 Performance Standard 21 CFR 1040.10 and 1040.11," as applicable.

#### Canadian EMC Standard

This instrument has been tested to and complies with ICES-001, Issue 4: "Industrial, Scientific, and Medical Radio Frequency Generators."

### European Safety, RoHS Directive, and EMC Standards

#### Safety

This instrument meets European requirements for safety (Low Voltage Directive 2014/35/EU). This instrument has been tested to and complies with standards EN61010-1, "Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use, Part 1: General Requirements."

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

EN 61010-2-081, "Particular Requirements for Automatic and Semi-Automatic Laboratory Equipment for Analysis and Other Purposes."

EN 60825-1, "Radiation Safety of Laser Products, Equipment Classification, Requirements, and User's Guide.

#### **EMC**

This instrument meets European requirements for emission and immunity (EMC Directive 2014/30/EU). This instrument has been tested to and complies with standard EN 61326 (Group 1, Class B), "Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements."

This instrument meets European RoHS Directive 2011/65/EU.

# Australian EMC Standards



This instrument has been tested to and complies with standard AS/NZS CISPR 11, "Limits and Methods Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radio-frequency Equipment."

Site Preparation Tasks

1

This chapter describes the tasks to perform before you install the Applied Biosystems StepOne<sup>TM</sup>or StepOnePlus<sup>TM</sup> Real-Time PCR System.

This chapter includes the following topics:

Overview	. 1-2
Assigning Personnel	1-3
Selecting the Site	. 1-4
Space Requirements	. 1-4
Environmental Requirements	1-7
Ventilation Requirements	. 1-8
Electrical Requirements	1-8
Computer Requirements	. 1-9
Network Requirements	1-10
Safety and Materials	1-11
Receiving and Inspecting the System	1-13
Moving the Crated Instrument to the Laboratory	1-14
During Installation	1-14

### Overview

Prepare your site for the installation according to the instructions in this chapter. Checklists are provided in Chapter 2, "Checklists."

# Site Preparation Schedule

To minimize the time between the shipment arrival and system installation:

- 1. Complete the site preparation tasks (Chapter 1).
- 2. Fill out the corresponding checklists (Chapter 2).

# Site Preparation Process

The general site preparation tasks and a suggested sequence for completing the tasks are summarized in Figure 1-1. The sequence can vary, but always:

- Review this guide first.
- Unpack and store the chemical installation kit as soon as you receive it.

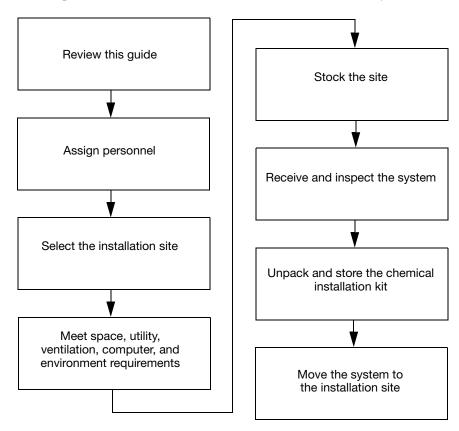


Figure 1-1 Site preparation tasks and their suggested sequence

# **Assigning Personnel**

### Laboratory Safety Representative

The laboratory safety representative should be familiar with laboratory safety procedures and know the location of all the safety equipment.

#### Tasks and Personnel

Table 1-1 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 1-1 Suggested personnel tasks

Personnel	Tasks
Site Preparation/ Installation Coordinator	<ul> <li>Reviews the site preparation guide for safety information and system requirements.</li> <li>Coordinates personnel and tasks.</li> <li>Orders required materials.</li> <li>Chooses the site.</li> <li>Reviews checklists with applicable personnel to verify that the site is properly prepared.</li> <li>Receives and inspects the system.</li> <li>Stores the Chemical Installation Kit.</li> <li>Schedules the installation and informs personnel of the installation date, if applicable.</li> <li>Ensures that the site is clear of unnecessary material on the installation day.</li> </ul>
Laboratory Safety Representative	<ul> <li>Reviews the site preparation guide for safety information.</li> <li>Ensures that the required safety practices and equipment are in place.</li> <li>Is available to assist with unpacking and setup.</li> </ul>
Laboratory Personnel/ Primary Users	<ul> <li>Review safety information.</li> <li>Ensure that all customer-provided materials for installation are present at the site.</li> </ul>
Facilities Personnel	<ul> <li>Ensure that installation requirements are met for:         <ul> <li>Space at the installation site</li> <li>Building clearances</li> <li>Temperature and humidity</li> <li>Electrical supply</li> <li>Computer</li> <li>Safety and installation materials</li> </ul> </li> <li>If possible, move the crated system to the site before the installation date.</li> </ul>
Network or IT Specialist (if the system will be connected to a network)	<ul> <li>Ensures that one active, tested local area network (LAN) connection is in place before the scheduled installation date.</li> <li>Ensures that network hardware is compatible with an RJ45-type connector.</li> <li>If necessary, supplies additional cables.</li> <li>Is available during installation to connect the system to the network.</li> <li>If applicable, provides and installs a network or dedicated printer.</li> </ul>

# Selecting the Site

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

- "Space Requirements" on page 1-4
- "Environmental Requirements" on page 1-7
- "Ventilation Requirements" on page 1-8
- "Electrical Requirements" on page 1-8
- "Computer Requirements" on page 1-9
- "Safety and Materials" on page 1-11

**IMPORTANT!** The site cannot be designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4) to be eligible for support. We do not install, service, or repair instruments in areas designated BSL-3 or BSL-4.

# **Space Requirements**

#### Instrument

Space requirements depends on your StepOne $^{TM}$  or StepOnePlus $^{TM}$  system purchase. Possible layouts:

- Standalone Instrument is not directly connected to a computer running StepOne<sup>™</sup> software. Use the standalone layout when the computer and the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument will be placed in separate locations.
- Colocated Instrument is connected to a computer by the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system cable. Use the colocated layout when both the computer and the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument will be placed together in the same location.

**Note:** You can choose to connect the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system to a network. See the *Applied Biosystems StepOne*<sup>TM</sup> and *StepOnePlus*<sup>TM</sup> *Real-Time PCR Systems Installation, Maintenance, and Networking Guide* on connecting the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system to an Ethernet network.



Figure 1-2 StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System

# System Components

If your purchase included a computer, then the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system includes the:

- StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument
- (Optional) Tower computer, monitor, keyboard, mouse, and control pad or

Laptop computer with mouse

**Note:** In some countries, the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system may be purchased without a computer. See your local representative for more information.

# Layout Requirements

#### Requirements:

- Avoid placing the system adjacent to heaters, cooling ducts, or in direct sunlight.
- Place away from any vibrating equipment such as a refrigerator or centrifuge. It should not share bench space with vibrating equipment.
- Place the computer within 2 m (6 ft) of the instrument in the colocated layout.
- Position the monitor, keyboard, and accessories to allow for proper ergonomics during use.

Figures 1-3 shows layout examples for the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System. For details on the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System space requirements, see Figure 1-4 on page 1-6 and Figure 1-5 on page 1-7.

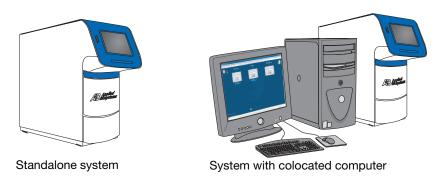


Figure 1-3 Examples of a standalone or colocated layout (not to scale)

# Dimensions and Weights

The table below indicates dimensions and weights of the system components, and computer components, if you ordered a computer system. Ensure that the installation site (floor space and/or bench space) can accommodate the dimensions and support the weights.

Component	Width	Depth	Height	Weight
StepOne <sup>TM</sup> or StepOnePlus <sup>TM</sup> system	24.6 cm (9.70 in.)	48.5 cm (19.11 in.)	51.80 cm (20.40 in.)	≈23.58±0.45 kg (53 ±1 lbs)
Computer, mini tower	18.7 cm (7.35 in.)	43.8 cm (17.24 in.)	41.1 cm (16.18 in.)	14.29 kg (31.50 lbs)

Component	Width	Depth	Height	Weight
Monitor	37.2 cm (14.64 in.)	18.2cm (7.16 in.)	45.8 cm (18.03 in.)	6.39 kg (14.08 lbs)
Keyboard	45.7 cm (18 in.)	17.8cm (7 in.)	5.1 cm (2 in.)	.68 kg (1.5 lbs)
Computer, laptop	33.8 cm (13.3_in.)	27.3 cm (10.75 in.)	3.58 cm (1.4 in.)	≈2.38 to 2.6 kg (5.24 to 5.73) lbs)

# Required Clearances

Required clearances for the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System are summarized below and illustrated in Figure 1-4 and Figure 1-5.

- Clearance on all sides At least 15.2 cm (6 in.) of clearance for ventilation, service access, and cable routing. Allow space for easy access to the back and sides.
- **Vertical clearance** At least 30.5 cm (12 in.) of unobstructed vertical clearance above the top of the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system to allow the top to be lifted during service.

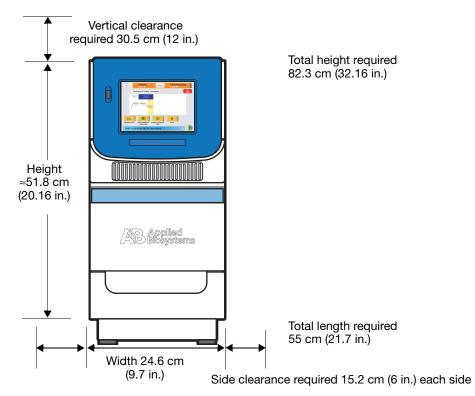


Figure 1-4 Space requirements front view (not to scale)

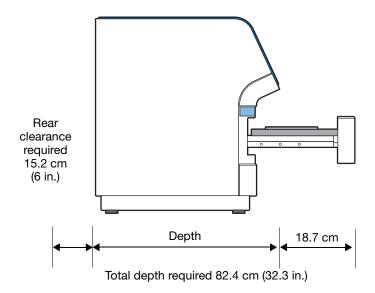


Figure 1-5 Space requirements side view (not to scale)

# **Environmental Requirements**

#### **Altitude**

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system is for indoor use only and for altitudes not exceeding 2000 m (6562 ft) above sea level.

# Temperature and Humidity Requirements

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

Condition	Acceptable Range
Temperature	15 to 30 °C (59 to 86 °F)  Maximum change of less than 15 °C per 24 hours
Temperature, transportation and storage	-30 to 60 °C (-22 to 140 °F)
Humidity	15 to 80% relative humidity, noncondensing
Humidity, transportation and storage	20 to 80% relative humidity

Fluctuations between day and night temperatures can cause system instability. For a stable environment, see "Laptop computer with mouse" on page 1-5.

#### **Pollution**

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System has a Pollution Degree rating of II. 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 II rating are laboratories and sales and commercial areas.

## **Ventilation Requirements**

### Venting Hot-Air-Only Exhaust

Hot-air exhaust is vented from the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system through the hot-air waste port on the back panel. The hot air exhaust is designed to dissipate heat produced by the instrument. The maximum thermal output of the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument is 1540 Btu/h (451.3 W). Consult your facilities department to determine if the laboratory ventilation system can maintain room temperature with this level of thermal output. If it can maintain room temperature during instrument operation, the hot-air exhaust port can be vented directly to room air.

# **Electrical Requirements**

#### Disconnecting Power

In case of emergency, you must be able to immediately disconnect the main power supply to the instrument.

# Power Connectors and Receptacles

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system is shipped to customers with 5 different power connectors to fit various country configurations. These connectors require standard 15 A wall receptacles with proper grounding. Do not use extension cords.

### Current Leakage Specification

The instrument meets the specification of 3- 3.5mA. Sites that test to ensure that current leakage is no more than 500 uA may want to purchase an isolation transformer to ground the instrument.

### System Electrical Requirements

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system can be configured for operating voltages between nominal voltages of 100VAC and 240VAC at 50 or 60 Hz. The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system is rated 900 Watts max. During installation, your installation coordinator configures the system for the proper input voltage.

**CAUTION**—Do not unpack or plug in any components until the appropriate person has configured the system for the proper operating voltage.

Table 1-2 provides electrical specifications for the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system. For all indicated input voltages, a 15 A circuit is required.

Table 1-2 Electrical specifications

Input Voltage (VAC)	Frequency (Hz)	Nominal Current Draw (A)	Power (W)
100	50/60	4.79	479
120	50/60	3.94	473

Table 1-2 Electrical specifications

Input Voltage (VAC)	Frequency (Hz)	Nominal Current Draw (A)	Power (W)
220	50/60	2.07	455
240	50/60	1.84	442

# Power Line Regulator

In areas where the supplied power is subject to voltage fluctuations exceeding  $\pm 10\%$  of the nominal value, a power line regulator may be required. High or low voltages can adversely affect the electronic components of the instrument.

# **Computer Requirements**

#### Minimum

We recommend the following minimum requirements for your computer system:

- Pentium IV processor
- USB port (1.1 or 2.0)
- 500 MHz
- 20 GB available on hard drive
- 512 MB of RAM
- Windows® XP operating system
- Computer:
  - UL 60950-1 listed laptop computer
- UL 60950-1 listed combination of a tower computer and monitor
- Ethernet network interface adaptor (10BASE-T)

# Software for Colocated Computer

The computer, if provided by Thermo Fisher Scientific, has the Microsoft<sup>®</sup> Windows<sup>®</sup> XP operating system with Service Pack 2 and Internet Explorer 6.x+ or Netscape Navigator 6.x+.

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 colocated computer additional software other than antivirus software. Changes to the configured software could void the instrument warranty and cause the system to be nonoperational.

# **Network Requirements**

#### **LAN Connection**

If the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System will be connected to a LAN, an active, tested LAN connection must be in place before your installation date. Due to differences in network connections, we cannot configure the system to access a specific network. For network issues, see your company IT department.

#### **Network Cables**

The computer if supplied is factory configured for the TCP/IP protocol. The product includes a fast Ethernet adapter (10/100baseT) with an RJ45-type connector and one 1.8 -m (6 -foot) Ethernet cable.

# Printer Requirements

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System can use a network or a dedicated printer. A dedicated printer is directly connected to the colocated computer.

# Safety and Materials

### **Safety Practices**

**IMPORTANT!** The site must not be designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4) to be eligible for Thermo Fisher Scientific support. We do not install, service, or repair 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. We expect 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 installation personnel will work.
- Appropriate fire extinguisher:
  - You are responsible for providing an appropriate fire extinguisher for use on or near the 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 MSDSs.

#### Materials for Installation

Provide the following materials for the installation:

- Safety glasses, lab coats, chemical-resistant, disposable gloves (powder-free)
- Lint-free tissues
- Ethanol, HPLC-grade or better
- 10% Bleach solution
- Water, Milli-Q® grade
- Three sizes of micropipettors and tips
  - $-1- to 10-\mu L$
  - -10- to  $100-\mu L$
  - -100- to 1,000- $\mu$ L
- Mini vortexer, centrifuge, and sample tubes

### Materials for Routine Operation

StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system supplies and consumables are necessary for routine operation of the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system. Before the system is installed, contact the sales representative to order supplies.

# Receiving and Inspecting the System

# Shipped Contents

The following are included in all StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system shipments:

- StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument
- Accessories:
  - 48-well or 96-well AB plates
  - USB drive
- Chemical installation kit

The following are included if you ordered them:

- Computer, monitor, computer keyboard, mouse, and control pad or
- Computer laptop with mouse

# Shipped Documentation

Refer to "How to Obtain More Information" on page 6 for a listing of documentation.

#### **Shipping List**

Verify that the items shown on the shipping list are the same items that you ordered.

### Inspecting Shipping Containers for Damage

Carefully inspect the shipping containers if the shipment was mishandled during transit and report any damage to customer support. Record any damage or mishandling on the shipping documents.

### Unpacking and Storing the Chemical Installation Kit

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

WARNING CHEMICAL HAZARD. Some chemicals used with Thermo Fisher Scientific instruments are potentially hazardous and can cause injury, illness, or death. Read and understand the Material Safety Data Sheets (MSDSs) provided by the chemical manufacturer before you store, handle, work with, or dispose of any chemicals or hazardous materials.

# Moving the Crated Instrument to the Laboratory

#### **Moving Schedule**

Before the date of installation:

- Clear the installation site of all unnecessary materials.
- If possible, have two people move the crated StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument from the receiving area to the installation site.
- 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 Applied Biosystems StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> Real-Time PCR System shipment contains the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> 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	614 cm (24 3/16 in.)
Length	663 cm (26 1/16 in.)
Depth	608 cm (23 15/16 in.)

# Instrument Weight

The StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> instrument weighs approximately 23.58±0.45 kg (53±1 lbs).

### Moving and Lifting the Instrument

**CAUTION PHYSICAL INJURY HAZARD.** 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.

# **During Installation**

After your installation personnel uncrate the system, installation and testing of the StepOne<sup>TM</sup> or StepOnePlus<sup>TM</sup> system either by your designated installer takes approximately 4 hours.

This chapter includes the following topics:	
Overview	2-2
Personnel Checklist	2-2
Space and Layout Checklist	2-3
Environmental Checklist	2-3
Ventilation Checklist	2-3
Electrical Checklist	2-4
Computer and Networking Checklist	2-4
Safety Checklist	2-5
Materials Checklist	2-6
System Receipt and Inspection Checklist	2-7
Moving the Crated Instrument Checklist	2-7

### 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. In the following checklists, date each item after verifying its completion.

## **Personnel Checklist**

For more information, see "Assigning Personnel" on page 1-3.

Date Verified	Designated Personnel
	Site Preparation/Installation coordinator
	Laboratory safety representative
	Laboratory personnel:
	To ensure that customer-supplied materials are on hand
	Primary users to be trained during installation and to subsequently train other users
	Facilities personnel:
	To provide environmental, electrical, and computer site-preparation requirements
	Two people to 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 1-4.

Date Verified	Requirements
	Location is away from:
	Heating or cooling ducts.
	Vibrating equipment, such as a refrigerator or centrifuge on the same benchtop.
	Direct sunlight.
	Computer workspace, if applicable, allows for proper ergonomics during use.
	Location accommodates the dimensions and weights specified in "Dimensions and Weights" on page 1-5.
	Location meets the requirements specified in "Required Clearances" on page 1-6.

# **Environmental Checklist**

For more information, see "Environmental Requirements" on page 1-7.

Date Verified	Requirement
	The altitude does not exceed 2000 m (6500 ft) above sea level.
	The conditions specified in "Temperature and Humidity Requirements" on page 1-7 have been met.
	Pollution Degree II - Only nonconductive pollutants, if any, are present.

# **Ventilation Checklist**

For more information, see "Ventilation Requirements" on page 1-8.

Date Verified	Requirement
	Instrument Hot-Air Exhaust Venting
	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 system (see "Venting Hot-Air-Only Exhaust" on page 1-8) is vented directly into the room air.  A suitable venting device such as a fume hood or fume dust is available to vent the hot air exhaust from the instrument space.

# **Electrical Checklist**

For more information, see "Electrical Requirements" on page 1-8.

Date Verified	Requirement	
	The main power supply to the instrument can be immediately disconnected.	
	Appropriate grounded power receptacles are available.	

# **Computer and Networking Checklist**

For more information, see "Computer Requirements" on page 1-9.

Date Verified	Requirement		
Antivirus Software			
	Appropriate antivirus software is available for loading on the system computer.		
Networking			
(If the system will be connected to a network)			
	One active, tested LAN connection is available.		
Network hardware is compatible with an RJ45-type connector. See "Network Requirements" on page 1-10.			
	For a colocated computer - If the network connection is more than 1.8 m (6 ft) from the system, a Category 5 RJ45 cable of the required length is available.		
Printer			
	A network printer or a dedicated printer and necessary print drivers are available.		

# **Safety Checklist**

For more information, see "Safety Practices" on page 1-11.

Date Verified	Requirement  The site is not designated BioSafety level 3 (BSL-3) or BioSafety level 4 (BSL-4) to be eligible for installation, service, and repair.				
	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				
	MSDSs readily available				

# **Materials Checklist**

For more information, see "Safety and Materials" on page 1-11.

Date Verified	Requirement			
	Materials for General Installation			
	Safety glasses and lab coats			
	Chemical-resistant disposable gloves (powder free)			
	Lint-free tissues			
	Ethanol, HPLC-grade or better			
	10% Bleach solution			
Water, Milli-Q grade				
	Three sizes of micropipettors and tips:			
	• 1- to 10-µL			
	• 10- to 100-μL			
	• 100- to 1000-μL			
	A mini vortexer, centrifuge, and sample tubes			
Materials for Routine Operation				
	Materials for routine operation after the installation are available or have been ordered (see "Materials for Routine Operation" on page 1-12).			

# System Receipt and Inspection Checklist

For more information, see "Receiving and Inspecting the System" on page 1-13.

Date Verified	Action		
	Verified that items on the packing list are those that were ordered. Otherwise, reported to customer support discrepancies in the packing list.		
	Opened and stored the Chemical Installation Kit components as specified in the kit operating instructions.		
	Received the system and inspected the shipping containers for mishandling or damage.		
	Reported to customer support:		
	Any damage to the shipping containers		
	Tip indicators or shock indicators that show evidence of mishandling during transit		

# Moving the Crated Instrument Checklist

For more information, see "Moving the Crated Instrument to the Laboratory" on page 1-14.

Date Verified	Item		
	The measured building clearances can accommodate the StepOne <sup>TM</sup> or StepOnePlus <sup>TM</sup> system crate dimensions (see "Required Building Clearances" on page 1-14). If the crate dimensions exceed building clearances, contact customer support.		
ir Z li t	Moved all the <i>crated</i> equipment to the laboratory before the date of the 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. No attempt should be made to lift the instrument.		
	Cleared the installation site of all unnecessary materials.		

# Index

A	dimensions 1-14
antivirus software 1-10	reporting shipping damage or tipping 1-13
Applied Biosystems	customer feedback, on Applied Biosystems documents 7
contacting 7 customer feedback on documentation 7	D
Technical Support 7	damage, reporting 1-13
assumptions for using this guide 5	DANGER, description 8
n	dimensions of system components 1-5
В	documentation, related 6
biohazardous waste, handling 14	_
biosafety level 3 or 4 1-4	E
building clearances, required 1-14	electrical
_	checklist 2-4
C	requirements 1-8
CAUTION, description 8	safety 15 specifications 1-8
checklists	•
computer 2-4	electromagnetic compatibility standards. <i>See</i> EMC standards
electrical 2-4	EMC standards 17
environmental 2-3	environmental checklist 2-3
inspection 2-7 layout 2-3	environmental requirements 1-7
materials 2-6	ergonomics, safety 16
moving the instrument 2-7	2.80.10.11.00, 04.20, 10
personnel assignments 2-2	F
safety 2-5	•
site preparation 2-1 system receipt 2-7	facilities personnel, tasks 1-3
Chemical Installation Kit, unpacking and storing 1-13	fire extinguisher 1-11
chemical safety 12	0
chemical waste safety 14	G
clearances	General 15
building 1-14	guidelines
required 1-6	chemical safety 12
computer checklists 2-4	chemical waste disposal 14 chemical waste safety 14
computer network. See network requirements	Chemical waste safety 14
computer software 1-10	Н
consumables, ordering 1-12	
conventions	hazard icons. See safety symbols, on instruments
IMPORTANTS! 6	hazard symbols. See safety symbols, on instruments
in this guide 5 safety 8	hazards. See safety
conventions used in this guide 5	Help system, accessing 7
See also moving the crated system	
and morning me traced by bream	

	requirements
inspection checklist 2-7	antivirus software 1-10
installation	component clearances and positioning 1-6
conditions for postponing 1-2	electrical 1-8
scheduling 1-2	environmental 1-7 installation 1-14
time required 1-14	layout 1-5
installation category 15	list of sections 1-4
installation coordinator, tasks 2-3	physical clearances 1-6
	space 1-4, 1-5, 1-6
instrument operation, safety 12	temperature and humidity 1-7
IT specialist, tasks 1-3	•
L	S
laboratory personnel, tasks 1-3	safety 1-11
	before operating the instrument 11
laboratory safety representative 1-3	biological hazards 15
layout checklist 2-3	checklist 2-5
layout requirements 1-5	chemical 12
lifting, safety 1-14	chemical waste 14 conventions 8
	electrical 15
M	ergonomic 16
materials	guidelines 12, 14
checklist 2-6	instrument operation 12
for routine operation 1-12	moving and lifting 1-14
moving and lifting, safety 11	moving and lifting instrument 11
	moving/lifting 11
moving the crated system 1-14	repetitive motion 16
moving the instrument checklist 2-7	standards 17
MSDSs	workstation 16
description 13	safety equipment, availability 1-11
obtaining 13	safety information
MSDSs, obtaining 7	lifting 1-14
N	safety labels, on instruments 10
IN	safety practices, required 1-11
network (IT) specialist, tasks 1-3	safety standards 17
network requirements, cable 2-4	safety symbols, on instruments 9
	shipping crate. See Crate
0	shipping damage, reporting 1-13
online Help. See Help system	shipping list, verifying 1-5, 1-13
overvoltage category (rating) 15	shipping, reporting damage 1-13
	site preparation
P	See also requirements
	checklists 2-1 flowchart 1-2
personnel assignments 1-3, 2-2	overview 1-2
personnel checklist 2-2	schedule 1-2
personnel tasks 1-3	selecting site 1-4
positioning system components 1-5	tasks 1-3
D	space checklist 2-3
R	space requirements 1-4, 1-5, 1-6
radioactive waste, handling 14	specifications, electrical 1-8
repetitive motion, safety 16	standards
reporting, shipping damage 1-13	EMC 17
	safety 17

```
supplies, ordering 1-12
symbols, safety 9
system
   building clearances 1-14
   components 1-4
   dimensions 1-5
   layout 1-5
   positioning components 1-5
   requirements 1-4
   weight 1-5
system receipt checklist 2-7
Т
tasks
   facilities personnel 1-3
   IT specialist 1-3
   laboratory personnel 1-3
   laboratory safety representative 1-3
   network specialist 1-3
   personnel 1-3
Technical Support, contacting 7
temperature and humidity, acceptable range 1-7
text conventions 5
training, information on 7
unpacking, Chemical Installation Kit 1-13
W
WARNING, description 8
waste disposal, guidelines 14
waste profiles, description 14
weight of system components 1-5
workstation safety 16
```

