



# Biacore™ X100

## Site Preparation Guide



# Table of Contents

<b>1</b>	<b>Introduction .....</b>	<b>3</b>
<b>2</b>	<b>Site preparation .....</b>	<b>5</b>
<b>3</b>	<b>Site preparation checklist .....</b>	<b>8</b>

# 1 Introduction

## Purpose of this document

The purpose of this document is to provide the information you need to fully prepare your site for the installation and testing of the system.

A site preparation checklist is included in this document.

## Scope of this document

This Site Preparation Guide covers the Biacore™ X100 instrument (product code BR110073).

## Installation by Cytiva

Read the entire Site Preparation Guide and complete the site preparation checklist in Chapter 3 before the product installation, to make sure that the site and personnel meet the installation requirements.

It is mandatory to return the completed checklist to Cytiva when booking an installation appointment.

## Contact Cytiva

To get information about returning the checklist, contact your Cytiva representative or follow the steps below to contact Cytiva.

Step	Action
1	Go to <a href="https://cytiva.com/contact">cytiva.com/contact</a> .
2	Select the applicable country from the <b>Change country</b> drop-down list.  <b>Note:</b> <i>If the applicable country is not listed, select the nearest region.</i>
3	In <b>Step 1</b> , select <b>I need help with an existing order</b> and click <b>Next</b> .
4	In <b>Step 2</b> , select <b>Order Issue &amp; Order Support</b> and click <b>Next</b> .
5	In <b>Step 3</b> , click <b>Message here</b> .  <i>Result:</i> The window with the contact form opens.
6	Fill in the required fields and click <b>Send</b> .

Step	Action
	<p><i>Result:</i> Your inquiry is sent to the Customer service team.</p>

## Assign personnel for site preparation tasks

The site preparation tasks listed in the following table can be assigned to more than one person at the customer site.

It is recommended that the same people assist the Cytiva service personnel during the subsequent installation process.

Personnel	Site preparation tasks
Site preparation/installation coordinator	<ul style="list-style-type: none"> <li>• Review the site preparation guide for safety information and system requirements.</li> <li>• Choose the site.</li> <li>• Coordinate personnel and tasks.</li> <li>• Coordinate the installation timing of the system with any domestic safety procedures and testings.</li> <li>• Order the required materials.</li> <li>• Gather information about the completed tasks, fill in the site preparation checklist, and return it to Cytiva before installation.</li> <li>• Schedule the installation and inform the relevant personnel of the installation date.</li> </ul>
Laboratory personnel/ Primary users	<ul style="list-style-type: none"> <li>• Review the safety information.</li> <li>• Make sure that all customer-provided materials for the installation are present at the site.</li> </ul>
Facilities personnel	<ul style="list-style-type: none"> <li>• Make sure that the installation requirements are met for: <ul style="list-style-type: none"> <li>- Space at the installation site</li> <li>- Environmental conditions</li> <li>- Biosafety</li> <li>- Ventilation and waste collection</li> <li>- Electrical supply</li> <li>- Safety and installation materials</li> </ul> </li> </ul>
IT support	<ul style="list-style-type: none"> <li>• Make sure that the computer requirements are fulfilled if the system controller is not from Cytiva. Refer to the <i>Biacore X100 data file</i> for more information.</li> <li>• Provide local administrator rights for the service engineer to install software at the time of the instrument installation.</li> </ul>

## 2 Site preparation

### Introduction

Before the product is delivered, the site must be prepared according to the information in this chapter.

### Select the site

Select a site that complies with national and international requirements for biosafety.

### Environmental conditions

The following requirements must be fulfilled at the location where the Biacore X100 system will be installed:

- Biacore X100 is intended for indoor use only
- The system units must not be exposed to direct sunlight
- The system units must not be placed adjacent to heaters or air-conditioners

### Environmental requirements

The installation site must comply with the following specifications:

Parameter	Specification
Ambient temperature, operation	18°C to 33°C
Relative humidity, operation	20% to 80%, non-condensing
Altitude, operation	Up to 2000 m

### Space requirements

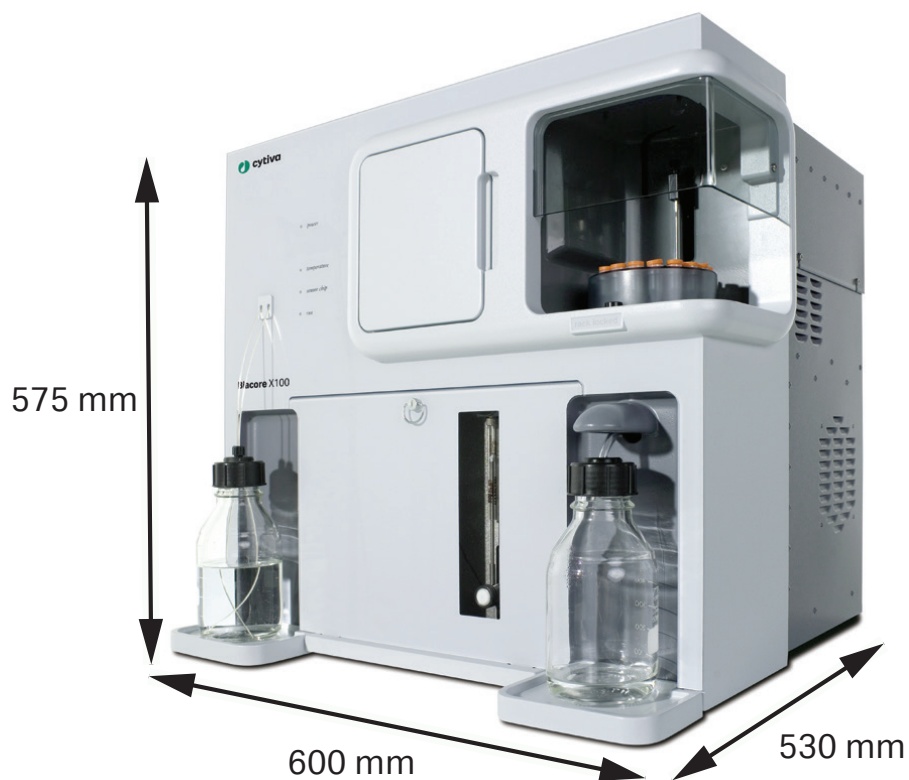
The following space requirements must be fulfilled at the location where the Biacore X100 system will be installed:

Parameter	Specification
Minimum bench area for operating the Biacore X100	100 x 75 cm

Parameter	Specification
Additional bench area for the computer with the control software	Approximately 50 x 75 cm
Clearance	At least 20 cm clearance is required on all sides of the instrument to allow adequate air circulation.

## Dimensions and weight

The following illustration shows the dimensions of the Biacore X100 instrument.



The Biacore X100 instrument weighs 47 kg.

## Electrical requirements

Parameter	Specification
Supply voltage	100 to 240 VAC
Frequency	50/60 Hz

Parameter	Specification
Maximum power	400 VA
Transient overvoltages	Overvoltage category II
Power outlet	Grounded
Number of wall outlets	3 Biacore X100 instrument, and system controller (computer and screen). See the manufacturer's specification for power requirements.

## Computer requirements

A computer supplied by Cytiva is recommended. If you do not purchase it from Cytiva, refer to the *Biacore X100 datafile* for information about computer requirements.

Refer to the manufacturer's documentation for specifications and environmental requirements of the computer.

Refer to the *Biacore X100 data file* for information about the included database.

## User-supplied accessories and consumables

The following equipment must be supplied by the user:

Parameter	Specification
Personal Protective Equipment (PPE) as required, (e.g., gloves, lab coats, etc.)	Used during installation, operation and maintenance
Water	Distilled and filtered water

## Other product specifications

Parameter	Specification
Noise	≤ 70 dB(A)

# 3 Site preparation checklist

## Introduction

Read the previous chapters in this site preparation guide before completing the checklist.

Complete the checklist in this chapter to make sure that all necessary preparations have been made for the installation of your Biacore X100.

You must send the completed checklist to a Cytiva office **before** Cytiva service personnel can install your Biacore X100. See details in Chapter 1.

## Checklist

Requirement	YES	NO	N/A	Comment/specify
<b>General</b>				
Has the Biacore X100 system been delivered? If yes, specify when the product was received (month/year)?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Will the Cytiva service personnel be required to take site EHS (Environment, Health and Safety) or security training before doing the installation?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Assigned personnel</b>				
Are personnel assigned as specified?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Will the personnel require training on how to use the Biacore X100 system?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Site selection</b>				
Does the site comply with biosafety requirements?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Are the general site requirements fulfilled?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Environmental requirements</b>				
Are the specified environmental conditions fulfilled?				
Ambient temperature	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Relative humidity	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	



Requirement	YES	NO	N/A	Comment/specify
Altitude	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Space requirements</b>				
Is the space available as specified?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Electrical requirements</b>				
Are the electrical requirements fulfilled as specified?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Number of wall outlets (3): <ul style="list-style-type: none"> <li>• One outlet for the Biacore X100 instrument</li> <li>• One outlet for the system controller computer</li> <li>• One outlet for the system controller screen</li> </ul>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>User supplied equipment</b>				
Personal Protective Equipment (PPE) as required, (e.g., gloves, lab coats, etc.).	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>User supplied consumables</b>				
Distilled and filtered water.	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	

### Additional information

### Customer contact information

<b>Date</b> (dd-mmm-yyyy)	<input type="text"/>		
<b>Company/Customer name</b>	<input type="text"/>		
Name	<input type="text"/>		
Address	<input type="text"/>		
Email	<input type="text"/>		
Phone number	<input type="text"/>		
<b>Customer experience level</b>	Beginner <input type="checkbox"/>	Intermediate <input type="checkbox"/>	Advanced <input type="checkbox"/>

Page intentionally left blank



cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

Biacore is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Any other third-party trademarks are the property of their respective owners.

© 2020–2023 Cytiva

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

29145041 AC V:7 05/2023