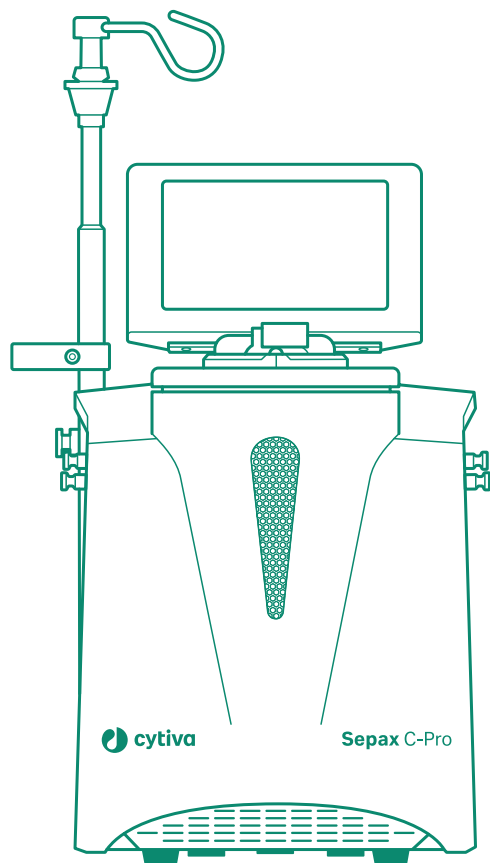


# Sepax™ C-Pro

## Site Preparation Guide



# Table of Contents

**1    Introduction ..... 3**

**2    Site preparation ..... 6**

**3    Site preparation checklist ..... 10**

# 1 Introduction

## Purpose of this document

The purpose of this document is to provide the information you need to prepare your site for the installation of the product.

A site preparation checklist is included in this document.

**Note:** *You must return the completed Site preparation checklist in Chapter 3 to Cytiva before Cytiva service personnel can visit your site to install your product.*

## Scope of this document

This Site Preparation Guide covers the Sepax™ C-Pro instrument.

## Installation by customer

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

## 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

Contact your Cytiva representative or follow the steps below to contact Cytiva about returning the checklist.

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> .

Step	Action
5	<p>In <b>Step 3</b>, click <b>Message here</b>.</p> <p><i>Result:</i> The window with the contact form opens.</p>
6	<p>Fill in the required fields and click <b>Send</b>.</p> <p><i>Result:</i> Your inquiry is sent to the Customer service team.</p>

## Assign personnel for site preparation tasks

The site preparation and installation tasks are typically split up between different persons and can take several weeks to prepare, depending on your organization and IT structure. It is up to the site preparation/installation coordinator to:

- Assign personnel and make sure that they are provided with sufficient information to complete the tasks, see table below.  
It is recommended that the same people assist the Cytiva service personnel during the subsequent installation process.
- Gather information about the completed tasks and fill in the site preparation checklist, see [Chapter 3 Site preparation checklist, on page 10](#).
- Return the completed checklist to Cytiva before installation, if applicable.

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>• Schedule the installation and inform the relevant personnel of the installation date.</li> <li>• Make sure the instrument box is not opened before the Cytiva service personnel arrive on site.</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> <li>- Lifting and moving equipment, if applicable</li> </ul> </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 general requirements must be fulfilled:

- The room must have exhaust ventilation.
- The instrument should not be exposed to sources of heat such as direct sunlight.
- Dust in the atmosphere should be kept to a minimum.

### Environmental requirements

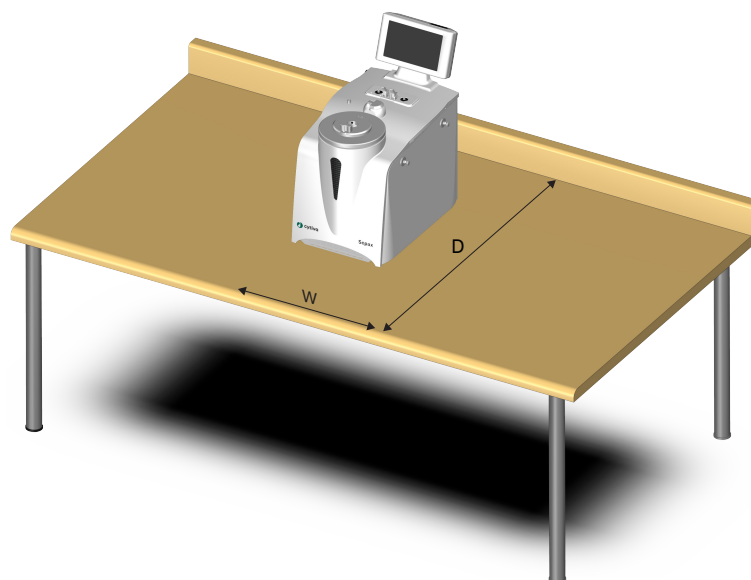
The installation site must comply with the following specifications.

Parameter	Requirement
Allowed location	Indoor use only
Ambient temperature, operating	15°C to 27°C
Ambient temperature, storage	0°C to 50°C
Relative humidity, operating	30% to 75%, non-condensing
Relative humidity, non-operating	20% to 75%, non-condensing
Altitude, operating	Up to 2000 m

### Space requirements

Sepax C-Pro instrument can be installed on a standard laboratory bench or on a trolley with wheels.

Prepare a clean working area on a stable laboratory bench or trolley that complies with the specifications in the following table. There must be sufficient space to turn the instrument on the bench for service access to the rear panel.



Parameter	Specification
Minimum bench area for operating the instrument (W x D)	47 x 70 cm
Free space required around the instrument	At least 20 cm free space in front of the instrument 10 cm free space on all other sides
Load capacity	980 N (100 kg) or higher
Inclination of bench surface	Horizontal $\pm 2^\circ$

## Dimensions and weight

The following illustration shows the dimensions of the instrument.



Parameter	Value
W	27 cm
D	40 cm
H1	46 cm
H2	96 cm
Weight	17 kg

## Electrical power requirements

The following table specifies the power requirements for the Sepax C-Pro instrument. For power requirements for the computer, refer to the manufacturer's specifications.

Parameter	Requirement
Supply voltage	100 to 240 V~ autorange
Frequency	50/60 Hz
Power outlet type	Grounded mains outlets
Transient overvoltages	Overvoltage category II



Parameter	Requirement
Power consumption	200 VA
Location of sockets	Maximum 2 m from the instrument (due to length of mains cable).

## Customer supplied equipment

The customer must supply the following:

- Personal Protective Equipment (PPE) as required.
- Lifting and moving equipment as required (refer to the *Operating Instructions*).
- Power plug for the mains cable.
- Equipotential bonding connection outside equipment to plant ground with a cross-sectional area of at least 4 mm<sup>2</sup>.

# 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 the product.

Send the completed checklist to a Cytiva office **before** Cytiva service personnel can install the Sepax C-Pro. See details in Chapter 1.

## Checklist

Requirement	YES	NO	N/A	Comment/specify
<b>General</b>				
Has the product been delivered)? If yes, specify when the product was received (month/year)?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Will 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>				
Have personnel been assigned as specified? See <a href="#">Assign personnel for site preparation tasks, on page 4</a>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Site selection</b>				
Does the site comply with requirements for biosafety?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Are the environmental conditions fulfilled?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Environmental requirements</b>				
Are the proper environmental conditions fulfilled as specified? See <a href="#">Environmental requirements, on page 6</a>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
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
Atmospheric pressure	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Space requirements</b>				
Is the specified space available? See <a href="#">Space requirements, on page 6</a>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Electrical requirements</b>				
Are the electrical requirements fulfilled as specified? See <a href="#">Electrical power requirements, on page 8</a>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>Network requirements</b>				
Will the product be connected to a network?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
Is there on-site IT support available if needed?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>User supplied equipment</b>				
Is Personal Protective Equipment (PPE) provided as required, (e.g., gloves, lab coats, etc.)? See <a href="#">Customer supplied equipment, on page 9</a>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	
<b>IQOQ (optional)</b>				
Will the IQOQ binder be delivered?	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	

Additional information

Customer contact information

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

Page intentionally left blank

**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://cytiva.com/techdocfeedback) or scan the QR code.



## cytiva.com

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

Sepax 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.

© 2023 Cytiva

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

29702829 AA V:2 02/2023