

BioCel Environmental- Control System

User Guide

Notices

© Agilent Technologies, Inc. 2008-2009

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

User Guide Part Number

G5500-90003

December/2007

Contact Information

Agilent Technologies Inc.
Automation Solutions
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
USA

Technical Support: 1.800.979.4811
or +1.408.345.8011
service.automation@agilent.com

Customer Service: 1.866.428.9811
or +1.408.345.8356
orders.automation@agilent.com

European Service: +44 12081443513
euroservice.automation@agilent.com

Documentation feedback:
documentation.automation@agilent.com

Web: <http://www.agilent.com>

Acknowledgements

Microsoft and Windows are registered trademarks of the Microsoft Corporation in the United States and other countries.

Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses


The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

If software is for use in the performance of a U.S. Government prime contract or sub-contract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14

(June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Safety Notices

 **A WARNING notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a WARNING notice until the indicated conditions are fully understood and met.**

A CAUTION notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a CAUTION notice until the indicated conditions are fully understood and met.

Letter to our Customers

Dear Customer,

The Agilent Technologies acquisition of Velocity11 resulted in the following changes:

- Creation of Agilent Technologies Automation Solutions, formerly Velocity11
- Renaming of some Velocity11 products
- New Customer Service and Technical Support contact information
- New website address for product information

Please make a note of the following changes as they impact this user guide.

Velocity11 product name changes

| Velocity11 product name | Changes to ... |
|--|-----------------------------|
| Access2 Automated Microplate Loader | Automated Centrifuge Loader |
| Element Automation System | BioCel 900 System |
| IWorks Device Driver Programming Interface | VWorks DCL Interface |
| PlatePierce Seal Piercing Station | Microplate Seal Piercer |
| VCode Barcode Print and Apply Station | Microplate Barcode Labeler |
| Velocity11 Robot | 3-Axis Robot |
| VHooks Integration Interface | VWorks Hooks Interface |
| VPrep Pipetting System | Vertical Pipetting Station |
| VSpin Microplate Centrifuge | Microplate Centrifuge |
| VStack Labware Stacker | Labware Stacker |

New contact information

Documentation feedback: documentation.automation@agilent.com

Technical Support: 1.800.979.4811 or +1.408.345.8011
service.automation@agilent.com

Customer Service: 1.866.428.9811 or +1.408.345.8356
orders.automation@agilent.com

European Service: +44 12081443513
euroservice.automation@agilent.com

Web: <http://www.agilent.com>

Contents

- Preface** v
 - Who should read this guide vi
 - What this guide covers vii
 - Accessing Velocity11 user guides viii

- Chapter 1. Environmental-Control System overview** 1
 - Environmental-Control System description 2
 - Nitrogen requirements 3
 - Hardware overview 4
 - Safety 7

- Chapter 2. Using the Environmental-Control System** 9
 - Workflow 10
 - Turning on and off the Environmental-Control System power 11
 - Setting the percentage of relative humidity 13
 - Closing and opening the doors 15
 - Turning on and off the nitrogen 17
 - Monitoring the percentage of relative humidity 19
 - Purging the nitrogen 20

- Appendix A. Schematic diagram** 21
 - Nitrogen and power system 22

Preface

This preface contains the following topics:

- “Who should read this guide” on page vi
- “What this guide covers” on page vii
- “Accessing Velocity11 user guides” on page viii

Who should read this guide

About this topic

This topic describes the target audience of this user guide.

Job roles

This user guide is for people with the following job roles:

| Job role | Responsibilities |
|---|--|
| Lab manager, administrator, or technician | Someone who is responsible for: <ul style="list-style-type: none"> <input type="checkbox"/> Managing the Environmental-Control System <input type="checkbox"/> Developing the applications that are run on it <input type="checkbox"/> Solving the more challenging problems that might arise <input type="checkbox"/> Developing training materials and standard operating procedures for operators |
| Operator | Someone who performs the daily production work on the Environmental-Control System and solves routine problems. Your organization may choose to create its own procedures for operators including the procedures in this guide. |

Related topics

| For more information about... | See... |
|--|---|
| What this guide covers | “What this guide covers” on page vii |
| How to access different formats of this user guide | “Accessing Velocity11 user guides” on page viii |

What this guide covers

What is covered This guide describes the Environmental-Control System in the BioCel[®] Automation System, the operation of the relevant hardware components, and the use of the iSeries Diagnostic software.

What is not covered This guide does not provide instructions for the BioCel System. For information about the BioCel System, see the *BioCel System User Guide*.

Software version This guide documents iSeries Diagnostics version 3.0.3 or later.

Related guides The *BioCel Environmental-Control System User Guide* should be used in conjunction with the following documents:

- BioCel User Guide*. The document explains how to use the system.
- Velocity11 automation control software user guides, such as the *VWorks Automation Control User Guide*. These user guides explain how to use the software to operate the system.

Related topics

| For more information about... | See... |
|--|---|
| Who should read this guide | “Who should read this guide” on page vi |
| How to access different formats of this user guide | “Accessing Velocity11 user guides” on page viii |

Accessing Velocity11 user guides

About this topic This topic describes the different formats of Velocity11 user information and explains how to access the user information.

Formats available Velocity11 user information is provided to you as:

- Online help
- A PDF file
- A printed book

The information in each format is the same but each format has different benefits.

Where to find user information

Online help

The online help is added to your computer with the Velocity11 lab automation system software installation.

PDF file

The PDF file of the user guide is on the software CD that is supplied with the product.

Velocity11 website

You can search the online help or download the latest version of any PDF file from the Velocity11 website at www.velocity11.com.

Note: All Velocity11 user information can be searched from the website at www.velocity11.com.

Online help

The online help is the best format to use when you are working at the computer and when you want to perform fast or advanced searches for information.

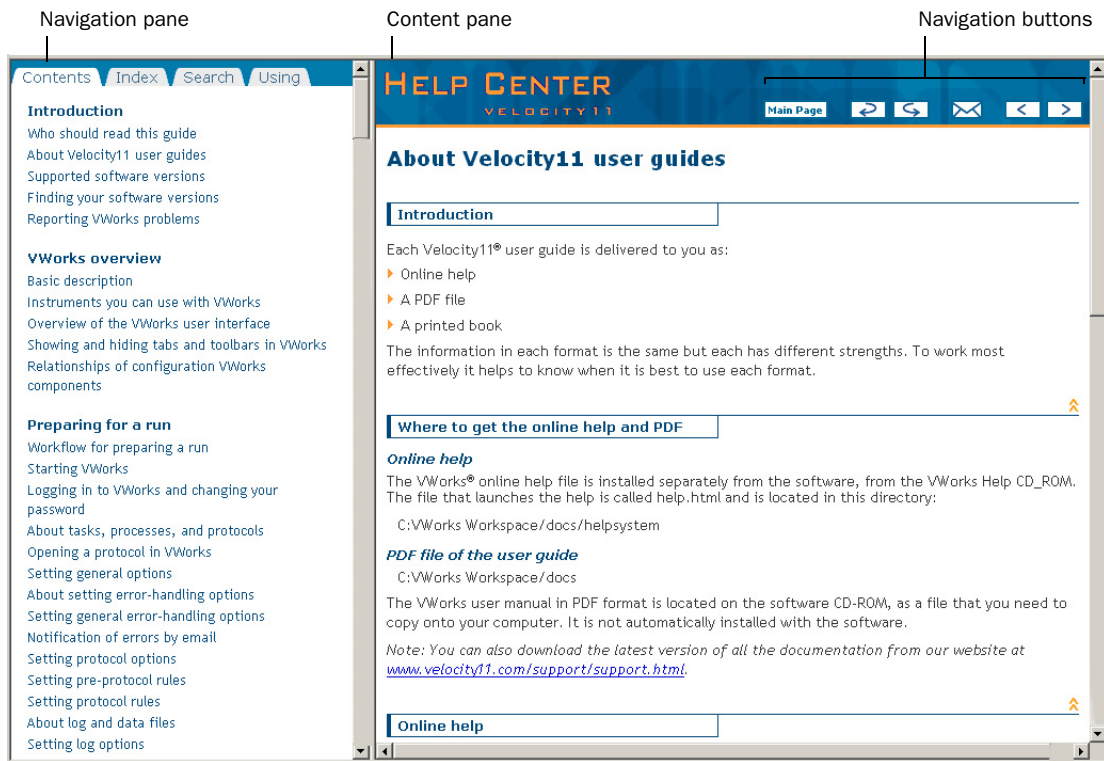
To open the online help:

In the Velocity11 lab automation software, press F1. The online help window opens.

Main features

The online help window contains the following:

- Navigation pane.* Consists of four tabs. The Contents, Index, and Search tabs provide different ways to locate information. The Using tab contains information about using the help system.
- Content pane.* Displays the online help topics.
- Navigation buttons.* Enables you to navigate through the pages. The online help includes a navigation pane, content pane, and navigation buttons.



PDF user guides

Computer requirements

To open a user guide in PDF format, you need a PDF viewer. You can download a free PDF viewer from the internet.

Printing and searching

The user guides in PDF format are mainly for printing additional copies. You can perform simple searches in the PDF file, although these searches are much slower than online help searches.

More information

For more information about using PDF documents, see the user documentation for the PDF viewer.

Related topics

| For more information about... | See... |
|-------------------------------|---|
| Who should read this guide | “Who should read this guide” on page vi |
| What this guide covers | “What this guide covers” on page vii |

Preface

BioCel Environmental-Control System User Guide

BioCel Environmental- Control System overview

1

This chapter contains the following topics:

- “Environmental-Control System description” on page 2
- “Nitrogen requirements” on page 3
- “Hardware overview” on page 4
- “Safety” on page 7

Environmental-Control System description

About this topic

This topic describes the BioCel Environmental-Control System and explains its use.

Description

The BioCel Environmental-Control System automatically maintains the desired percentage of relative humidity within the BioCel System. The system uses nitrogen gas to displace air inside the sealed BioCel System chamber, thereby reducing the humidity.

Components

The Environmental-Control System consists of the following components:

- Nitrogen On and Off buttons on the Environmental Controls panel
- Nitrogen Purge switch on the Environmental Controls panel
- Door-locking and interlock mechanisms
- Diagnostics software that allows you to set the desired percentage of relative humidity

Before you operate the system

!! INJURY HAZARD !! For safe operation, it is imperative that you follow the precautions in “Safety” on page 7.

Related topics

| For information about... | See... |
|--|-----------------------------------|
| Environmental-Control System specifications | “Nitrogen requirements” on page 3 |
| Environmental-Control System component description | “Hardware overview” on page 4 |
| Safety information | “Safety” on page 7 |

Nitrogen requirements

About this topic

This topic lists the nitrogen requirements for the Environmental-Control System. For the BioCel System physical, electrical, and environmental requirements, see the *BioCel System User Guide*.

Requirements

| Requirement | Value |
|-------------|--|
| Quality | Application dependent |
| Source | One of the following: <input type="checkbox"/> Centralized source (house) <input type="checkbox"/> Compressed-nitrogen cylinders |
| Pressure | 0.6 MPa (100 psig) maximum |
| Flow rate | 1100 Lpm (40 scfm) maximum |

Note: The flow rate you use depends on the time required to achieve the target percentage of relative humidity.

Connection to the BioCel System

Nitrogen lines are connected at the top of the BioCel System. For the location of the external utilities connections, see the *BioCel System User Guide*.

Related topics

| For information about... | See... |
|---|---------------------------------|
| Environmental-Control System components | “Hardware overview” on page 4 |
| Safety information | “Safety” on page 7 |
| BioCel System lab requirements | <i>BioCel System User Guide</i> |

Hardware overview

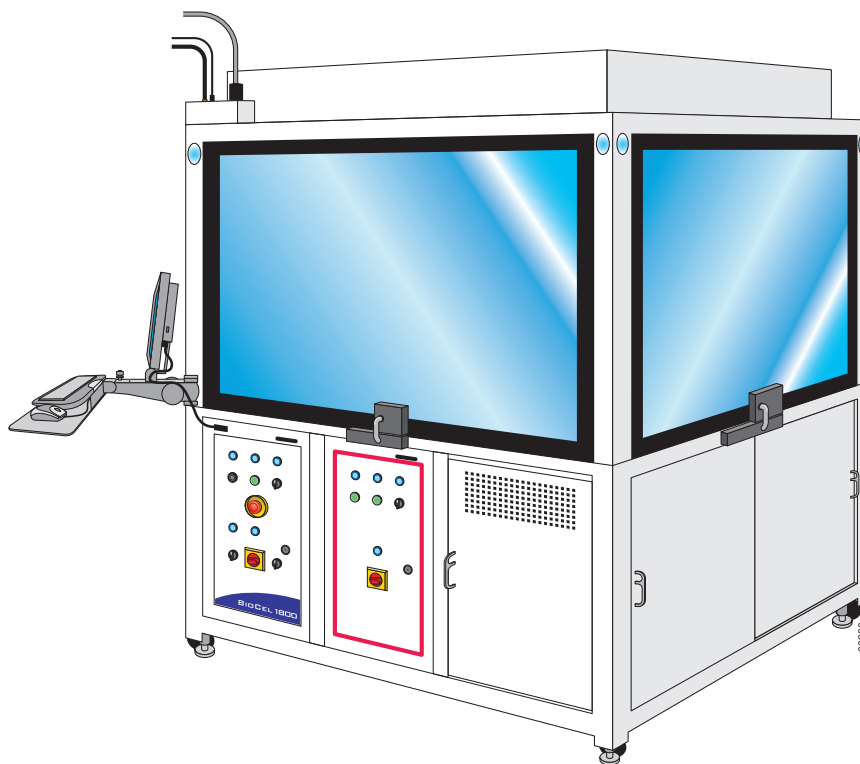
About this topic

This topic describes the following Environmental-Control System hardware features:

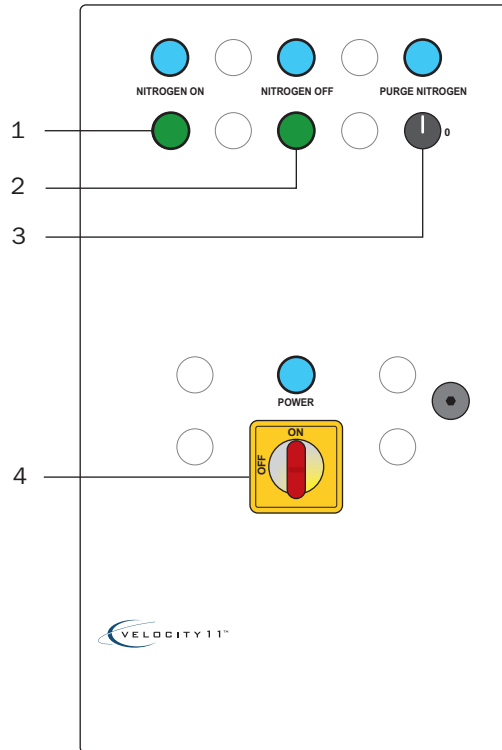
- Environmental Controls panel
- Nitrogen Utilities panel

Environmental Controls panel

The Environmental Controls panel consists of a switch for power, buttons for turning on and off the nitrogen, and a switch for purging the nitrogen. The Environmental Controls panel is located next to the power panel.



The following figure shows the Environmental Controls panel in detail.

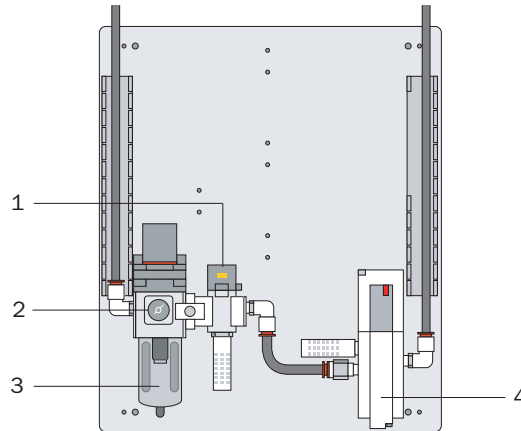


| Item | Feature | Description |
|------|-----------------------|--|
| 1 | NITROGEN ON button | Turns on the nitrogen to the system. |
| 2 | NITROGEN OFF button | Turns off the nitrogen to the system. |
| 3 | PURGE NITROGEN switch | Purges the nitrogen from the system chamber. When you turn the PURGE NITROGEN switch, the nitrogen is turned off automatically. |
| 4 | POWER switch | Turns on and off the Environmental-Control System. |

Nitrogen Utilities panel

The Nitrogen Utilities panel contains the nitrogen-supply connection, regulator, and filter. The utilities panel is located behind the Environmental Controls panel. To access the panel, slide open the cabinet door nearest the Environmental Controls panel.

The following figure shows the components that control the nitrogen supply in detail.



| Item | Feature | Description |
|------|---|---|
| 1 | Nitrogen supply shutoff valve and residual pressure release | Turns on or off the nitrogen supply to the system. Use the valve to turn off the nitrogen supply only if the NITROGEN ON or the NITROGEN OFF buttons on the Environmental Controls panel is not working, or if the system is being serviced. |
| 2 | Nitrogen pressure regulator | Controls the nitrogen pressure to the system. |
| 3 | Filter | Filters the air supply to the system. |
| 4 | Interlock air shutoff valve | Shuts off the air to moving parts when the interlock circuit is interrupted. |

Related topics

| For information about... | See... |
|------------------------------|--|
| Environmental Controls panel | "Environmental Controls panel" on page 4 |
| Nitrogen requirements | "Nitrogen requirements" on page 3 |
| Safety information | "Safety" on page 7 |

Safety

About this topic

!! INJURY HAZARD !! It is imperative that you follow the precautions in this section and in the *BioCel System User Guide* for safe operation.

The BioCel System with the Environmental-Control System is designed for safe operation. Under normal operating conditions, you are protected from high-pressure gas and high levels of nitrogen. However, you must be aware of the potential hazards and understand how to avoid being exposed to them.

This section presents the Environmental-Control System safety information. Read this section in conjunction with the safety information in the *BioCel System User Guide*.

Safety standards

The BioCel System with the Environmental-Control System is CE-certified and complies with the following CE safety directives: EN61010-1:1993, including Amendment 2:1995 (Safety).

High-pressure gas cylinders

Nitrogen gas is used to displace air inside the BioCel System.

Follow the local, state, and federal safety codes for the placement and mounting of gas cylinders. For example, you might have to attach a standard cylinder bracket to a solid permanent structure to meet or exceed all local seismic and safety requirements.

Always use good lab practices when handling high-pressure cylinders. Make sure you follow any instructions provided with the cylinders.

Nitrogen gas

!! INJURY HAZARD !! Nitrogen is an odorless, colorless, and nontoxic gas that can cause suffocation by displacing air.

Always use nitrogen in a well-ventilated area. Always turn off the nitrogen source when the system is not in use.

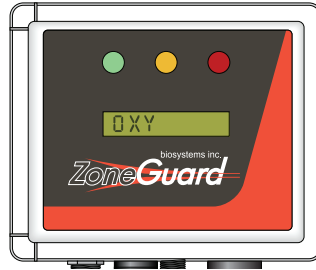
The following safety label is placed in the BioCel System to warn you of low-oxygen hazards.



Oxygen monitors

!! INJURY HAZARD !! Do not remove the provided oxygen monitors.

Two oxygen monitors are provided with the system: one on the inside of the system, and one on the outside of the system. The oxygen monitors sense and display the percentage of oxygen present.



To prevent accidental suffocation:

- The internal monitor interacts with the safety interlock system so that when the system chamber contains less than 19.5% oxygen, the doors lock automatically. To open the doors, the system requires that you first purge the nitrogen.
- The external monitor interacts with the safety interlock system so that when the air surrounding the BioCel System contains less than 19.5% oxygen, the shutoff valves cut off the nitrogen supply to the system.

Door locks and safety interlock system

!! INJURY HAZARD !! Do not disable or defeat the door locks or interlocks.

The BioCel System has a door-locking and safety interlock system that prevent you from opening the doors and being exposed to high nitrogen levels. When you turn on nitrogen at the Environmental Controls panel, the doors lock automatically. To unlock the doors, you must first purge the nitrogen.

Related topics

| For information about... | See... |
|------------------------------|--|
| Environmental Controls panel | "Environmental Controls panel" on page 4 |
| Nitrogen requirements | "Nitrogen requirements" on page 3 |
| Operating instructions | "Using the Environmental-Control System" on page 9 |

Using the Environmental-Control System

2

This chapter explains how to operate the Environmental-Control System. The topics are:

- “Workflow” on page 10
- “Setting the percentage of relative humidity” on page 13
- “Closing and opening the doors” on page 15
- “Turning on and off the nitrogen” on page 17
- “Purging the nitrogen” on page 20

Workflow

Workflow steps

The following table presents the steps for using the Environmental-Control System for a protocol run.

Note: Velocity11 sets up the profile for the Environmental-Control System. Do not change the existing profile settings.

| Step | For this task... | See... |
|------|--|---|
| 1 | Start up the BioCel System. | <i>BioCel System User Guide</i> |
| 2 | Turn on the Environmental-Control System. | “Turning on and off the Environmental-Control System power” on page 11 |
| 3 | Set the desired percentage of relative humidity. | “Setting the percentage of relative humidity” on page 13 |
| 4 | Close the doors. | “Closing and opening the doors” on page 15 |
| 5 | Turn on the nitrogen. | “Turning on and off the nitrogen” on page 17. |
| 6 | Start a protocol run. | Velocity11 automation control software user guide, such as the <i>VWorks User Guide</i> |
| 7 | Monitor the percentage of relative humidity. | “Setting the percentage of relative humidity” on page 13 |
| 8 | Purge the nitrogen. | “Purging the nitrogen” on page 20 |

Related topics

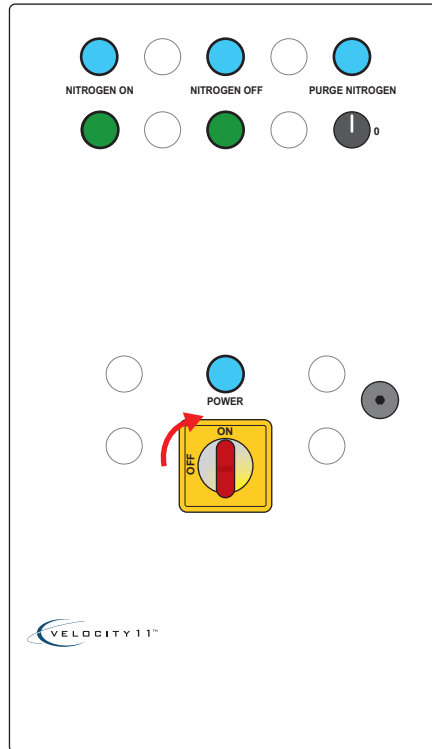
| For information about... | See... |
|---|---|
| Operating the BioCel System | <i>BioCel System User Guide</i> |
| Running a protocol | Velocity11 automation control software user guide, such as the <i>VWorks User Guide</i> |
| Environmental-Control System components | “Hardware overview” on page 4 |
| Nitrogen requirements | “Nitrogen requirements” on page 3 |
| Safety information | “Safety” on page 7 |

Turning on and off the Environmental-Control System power

Turning on the power

To turn on the Environmental-Control System:

On the Environmental Controls panel, turn the **POWER** switch clockwise.

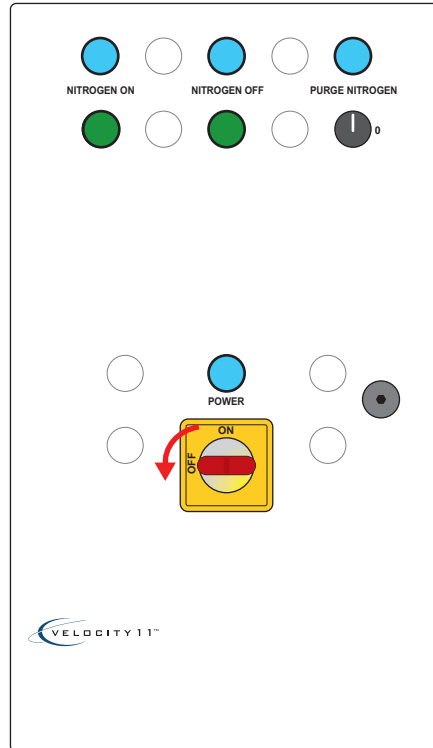


Turning off the power

After the protocol run is finished, turn off the nitrogen, and then turn off the Environmental-Control System.

To turn off the Environmental-Control System:

On the Environmental Controls panel, turn the **POWER** switch counterclockwise.



Related topics

| For information about... | See... |
|---|---|
| Set the desired percentage of relative humidity | "Setting the percentage of relative humidity" on page 13 |
| Closing or opening the doors | "Closing and opening the doors" on page 15 |
| Turning on or turning off the nitrogen | "Turning on and off the nitrogen" on page 17 |
| Monitoring the internal environment | "Monitoring the percentage of relative humidity" on page 19 |
| Purging the nitrogen | "Purging the nitrogen" on page 20 |
| Nitrogen requirements | "Nitrogen requirements" on page 3 |
| Safety information | "Safety" on page 7 |

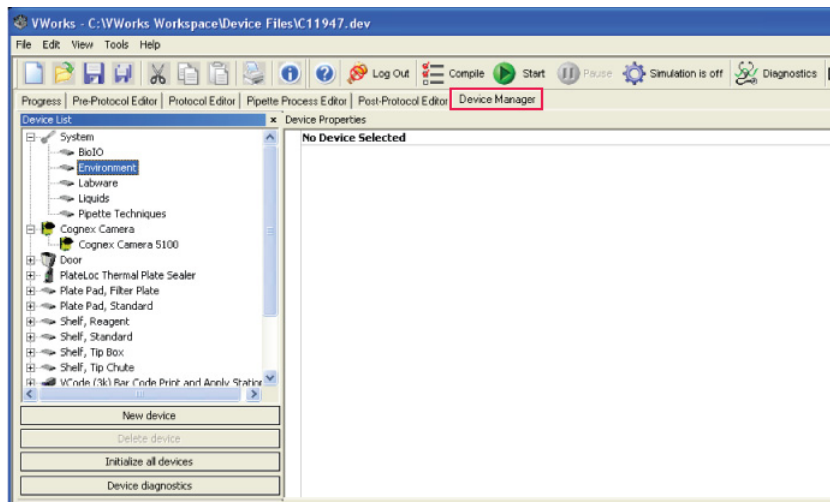
Setting the percentage of relative humidity

Procedure

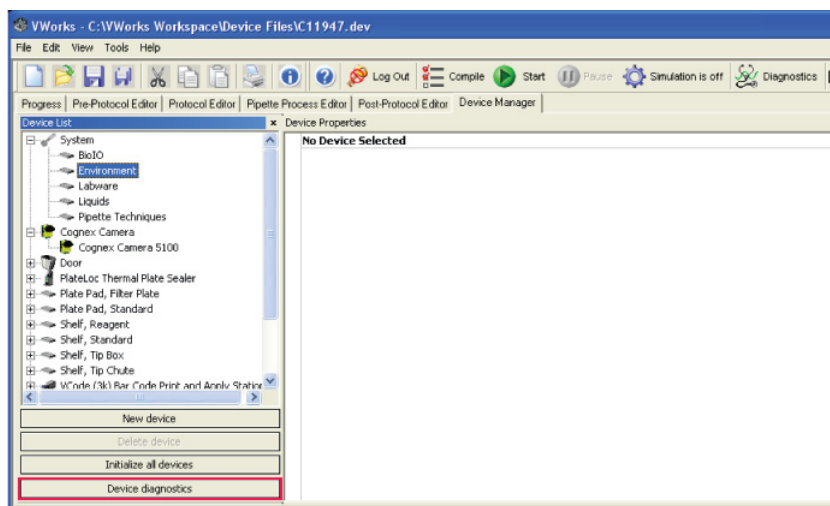
You set the desired percentage of relative humidity in the iSeries Diagnostics software.

To set the desired percentage of relative humidity:

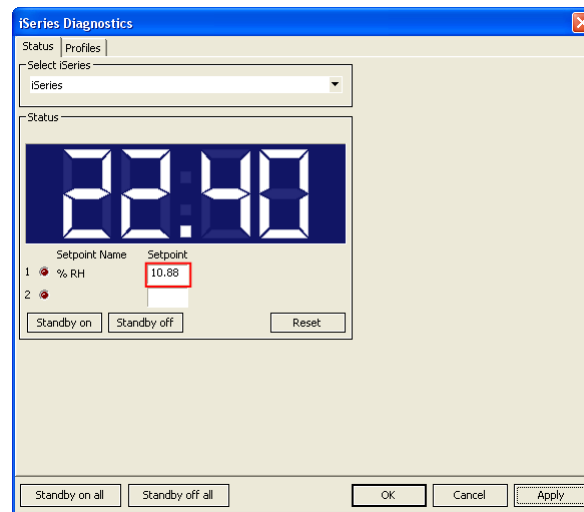
1. In the Velocity11 automation control software, click the **Device Manager** tab.



2. In the **Device List** area, select **Environment**, and then click **Device Diagnostics**. The iSeries Diagnostics dialog box opens.



- In the **Status** tab, type the desired percentage of relative humidity in the **% RH Setpoint** box.



- Click **OK** to save the setting and return to the Velocity11 automation control software window.

You can click **Apply** to apply the setting without saving it.

Related topics

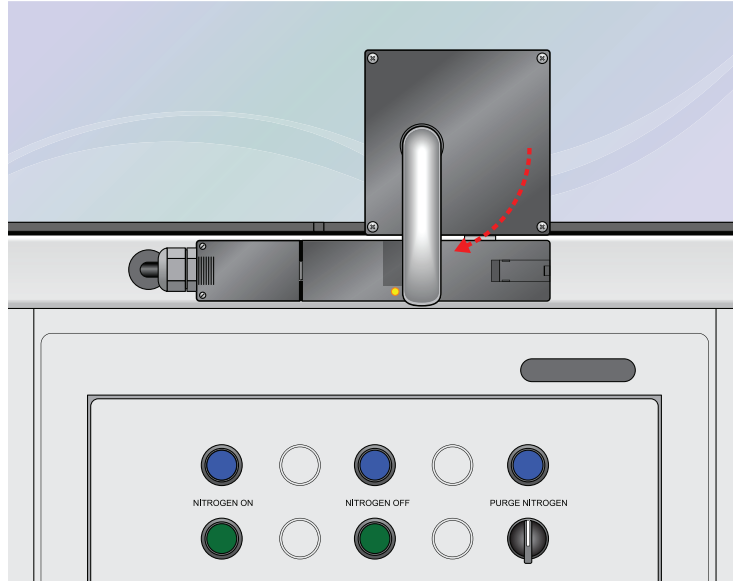
| For information about... | See... |
|---|--|
| Turning on the Environmental-Control System | “Turning on and off the Environmental-Control System power” on page 11 |
| Closing or opening the doors. | “Closing and opening the doors” on page 15 |
| Turning on or turning off the nitrogen | “Turning on and off the nitrogen” on page 17 |
| Nitrogen requirements | “Nitrogen requirements” on page 3 |
| Safety information | “Safety” on page 7 |

Closing and opening the doors

Closing the doors

To close the BioCel System doors:

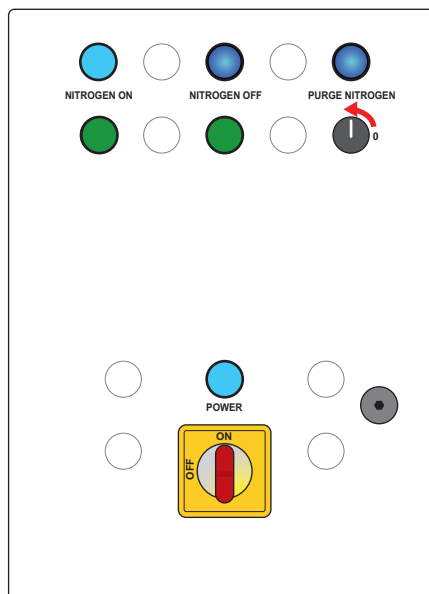
1. Close the system doors.
2. Turn the handle clockwise.



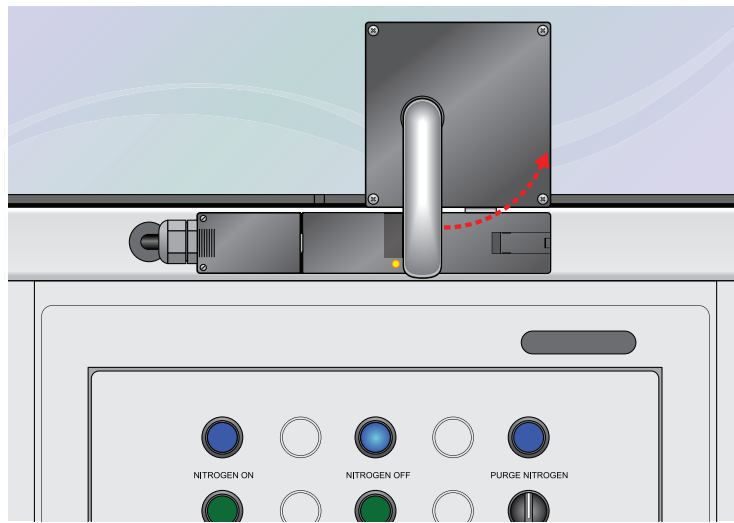
Opening the doors

To open the BioCel System doors:

1. On the Environmental Controls panel, turn the **PURGE NITROGEN** switch counterclockwise. The indicator light above the switch turns on. The nitrogen automatically turns off. The doors automatically unlock when the oxygen inside reaches a safe level.



- Turn the handle counterclockwise.



- Lift the door up.

Related topics

| For information about... | See... |
|---|--|
| Turning on the Environmental-Control System | “Turning on and off the Environmental-Control System power” on page 11 |
| Setting the desired percentage of relative humidity | “Setting the percentage of relative humidity” on page 13 |
| Turning on the nitrogen | “Turning on and off the nitrogen” on page 17 |
| Nitrogen requirements | “Nitrogen requirements” on page 3 |
| Purging the nitrogen | “Purging the nitrogen” on page 20 |
| Safety information | “Safety” on page 7 |

Turning on and off the nitrogen

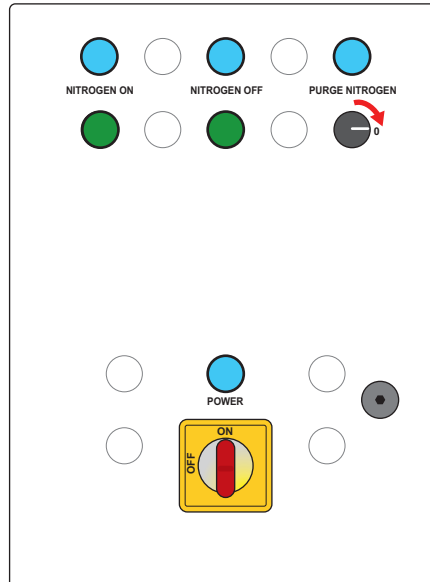
Before you start

Before you turn on the nitrogen, make sure the doors are closed and the handles are positioned vertically. See “Closing the doors” on page 15.

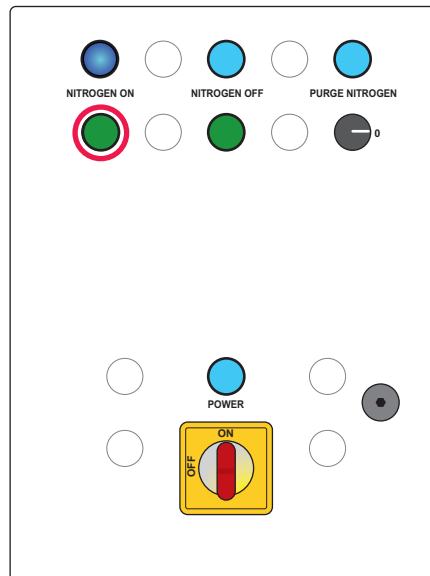
Turning on the nitrogen

To turn on the nitrogen:

1. Make sure the **PURGE NITROGEN** switch is in the off position (O). The indicator light above the switch should be off.



2. On the Environmental Controls panel, press the **NITROGEN ON** button. The indicator light above the button turns on.

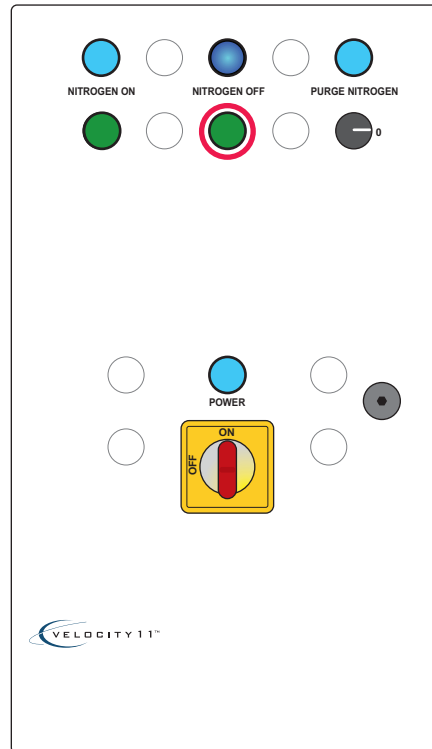


Turning off the nitrogen

The system turns off the nitrogen automatically when you press the PURGE NITROGEN button on the Environmental Controls panel.

To turn off the nitrogen manually:

On the Environmental Controls panel, press the **NITROGEN OFF** button. The indicator light above the button turns on.



Related topics

| For information about... | See... |
|---|--|
| Safety information | "Safety" on page 7 |
| Nitrogen requirements | "Nitrogen requirements" on page 3 |
| Turning on the Environmental-Control System | "Turning on and off the Environmental-Control System power" on page 11 |
| Setting the desired percentage of relative humidity | "Setting the percentage of relative humidity" on page 13 |
| Closing the doors | "Closing and opening the doors" on page 15 |
| Purging the nitrogen | "Purging the nitrogen" on page 20 |

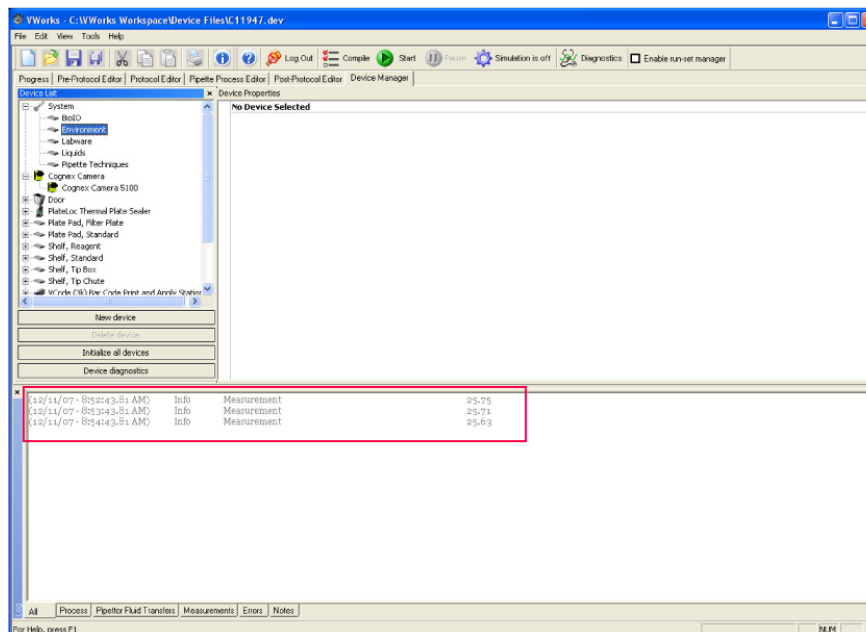
Monitoring the percentage of relative humidity

Procedure

The system checks and displays the percentage of relative humidity constantly.

To view the humidity level, in the Velocity11 automation control software window:

Check the measurements displayed in the Log area at the bottom of the window.



Related topics

| For information about... | See... |
|---|--|
| Turning on the Environmental-Control System | “Turning on and off the Environmental-Control System power” on page 11 |
| Setting the desired percentage of relative humidity | “Setting the percentage of relative humidity” on page 13 |
| Closing the doors | “Closing and opening the doors” on page 15 |
| Turning on the nitrogen | “Turning on and off the nitrogen” on page 17 |
| Nitrogen requirements | “Nitrogen requirements” on page 3 |
| Safety information | “Safety” on page 7 |

Purging the nitrogen

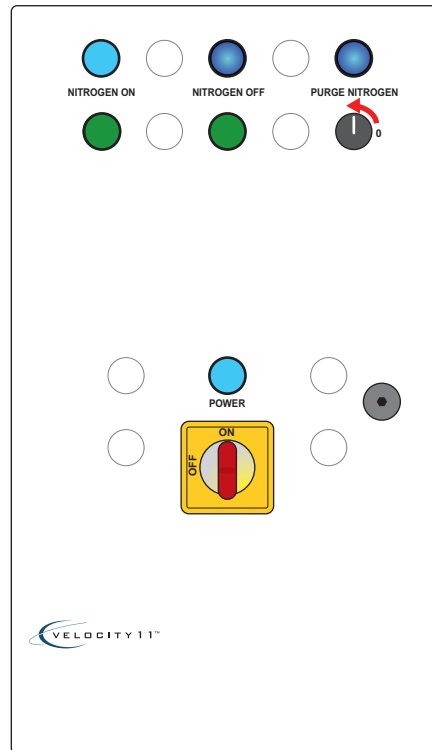
Procedure

The system locks the doors when the internal oxygen level is 19.5% or lower. To open the doors, you must first purge the nitrogen.

Note: The system purges the nitrogen automatically when the external oxygen monitor senses that the oxygen surrounding the BioCel System drops below 19.5%.

To purge the nitrogen manually:

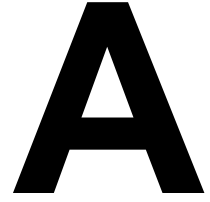
On the Environmental Controls panel, turn the **PURGE NITROGEN** switch counterclockwise. The indicator light above the switch turns on. The nitrogen turns off automatically. Fans in the system expel the nitrogen and draw air from outside. The system requires approximately 3 minutes for the internal oxygen level to reach 19.5%.



Related topics

| For information about... | See... |
|--------------------------|--|
| Opening the doors | “Closing and opening the doors” on page 15 |
| Turning off the nitrogen | “Turning on and off the nitrogen” on page 17 |
| Safety information | “Safety” on page 7 |

Schematic diagram



This appendix provides the schematic diagram and a description of the nitrogen system.

Nitrogen and power system

About this topic

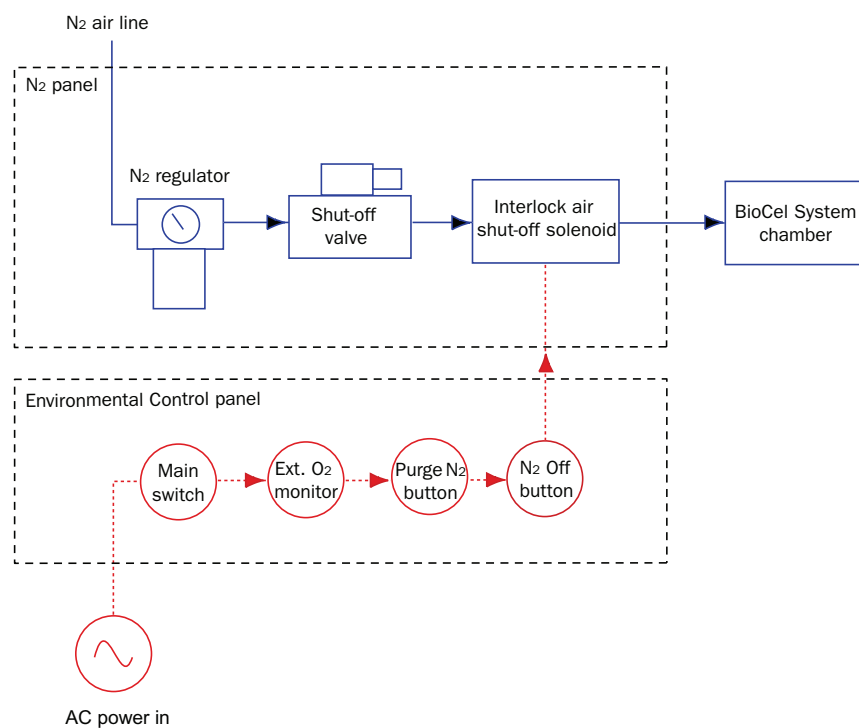
This topic summarizes the interaction between the nitrogen supply and power system. The information can be useful for troubleshooting purposes.

Diagram and description

The BioCel System uses nitrogen to displace air, thereby reducing the percentage of relative humidity. The flow of nitrogen through the system is controlled by the nitrogen pressure regulator in the nitrogen utilities panel (“Nitrogen Utilities panel” on page 6).

The following diagram shows when nitrogen is cut off from the system:

- The BioCel System is turned off (main switch).
- The external oxygen monitor senses the oxygen level dropped below 19.5%.
- The PURGE NITROGEN button is pressed.
- The NITROGEN OFF button is pressed.
- The nitrogen shutoff valve in the nitrogen utilities panel is turned off.



Related topics

| For information about... | See... |
|---------------------------------|--|
| Environmental Controls panel | “Environmental Controls panel” on page 4 |
| Nitrogen utilities panel | “Nitrogen Utilities panel” on page 6 |
| Power panel | <i>BioCel System User Guide</i> |
| Safety information | “Safety” on page 7 |



User Guide
G5500-90003