

ÄKTA go™

Operating Instructions

Original instructions

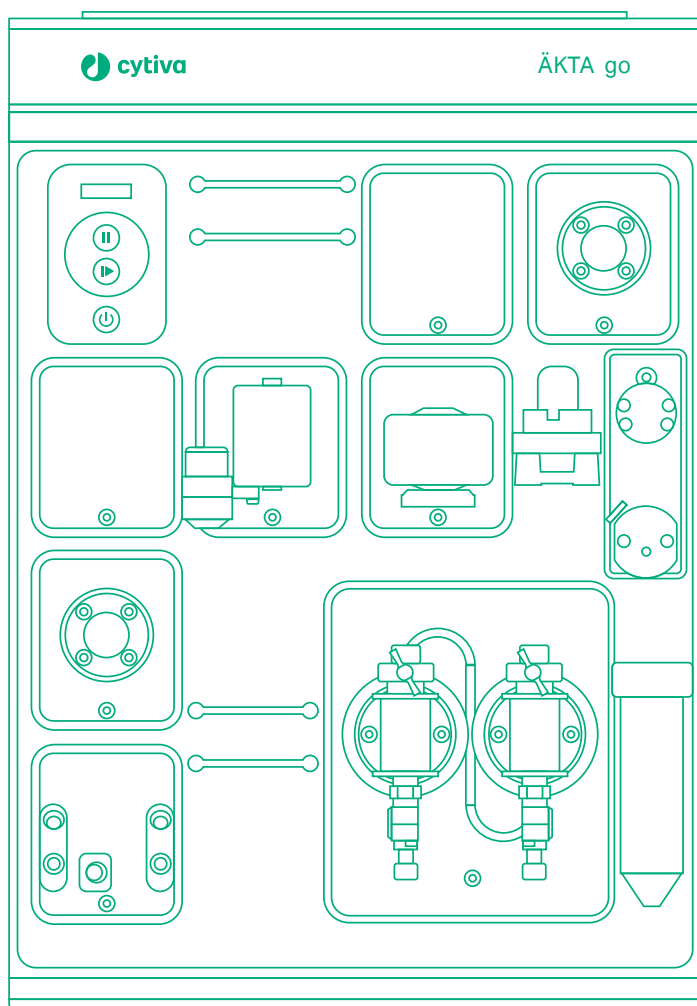


Table of Contents

1	Introduction	4
1.1	Important user information	5
1.2	About this manual	6
1.3	Associated documentation	7
1.4	Abbreviations	9
2	Safety instructions	10
2.1	Safety precautions	11
2.2	Labels and symbols	15
2.3	Emergency procedures	16
3	System description	18
3.1	ÄKTA go	19
3.2	Available modules	25
3.3	UNICORN	28
4	Installation	32
4.1	Site preparation	33
4.1.1	<i>Delivery, storage, and unpacking</i>	34
4.1.2	<i>Space requirements</i>	36
4.1.3	<i>Site environment</i>	39
4.1.4	<i>Power requirements</i>	41
4.1.5	<i>Computer requirements</i>	42
4.1.6	<i>Required materials</i>	43
4.2	Hardware installation	44
4.2.1	<i>Connect the system</i>	45
4.2.2	<i>Install waste tubing</i>	48
4.2.3	<i>Prepare the pump rinsing system</i>	50
4.2.4	<i>Start the instrument</i>	53
4.3	Software installation	54
4.3.1	<i>Download and Install UNICORN</i>	55
4.3.2	<i>Download the Instrument Configuration</i>	56
4.3.3	<i>Adapt UNICORN to your system</i>	57
4.4	Start UNICORN and connect to system	59
5	Prepare the system for a run	62
5.1	Prepare the flow path	64
5.2	Prime inlets and purge pump heads	66
5.3	Performance tests	72
5.4	Connect a column	75
5.5	Pressure alarms	78
5.6	Prepare for a run at low temperature	79
6	Run a method	80
6.1	Create a method	81
6.2	Prepare sample for loading	84
6.3	Start a method run	86

6.4	Monitor or interact with the run	87
6.5	Evaluate the run	88
6.6	Procedures after the run	90
7	Maintenance	93
7.1	Clean the instrument externally	94
7.2	Perform system Cleaning-In-Place (CIP)	95
7.3	Replace pump rinsing liquid	99
7.4	Replace the main fuses	100
8	Troubleshooting	102
9	Reference information	104
9.1	System specifications	105
9.2	Chemical resistance specifications	106
9.3	Recycling information	108
9.4	Regulatory information	109
9.4.1	Contact information	110
9.4.2	European Union and European Economic Area	111
9.4.3	Great Britain	112
9.4.4	Eurasian Economic Union (Евразийский экономический союз)	113
9.4.5	North America	115
9.4.6	General regulatory statements	116
9.4.7	South Korea	117
9.4.8	China	118
9.4.9	Health and Safety Declaration Form	121
Index		123

1 Introduction

This chapter contains information about this manual and associated user documentation, important user information and intended use of ÄKTA go™.

In this chapter

Section		See page
1.1	Important user information	5
1.2	About this manual	6
1.3	Associated documentation	7
1.4	Abbreviations	9

1.1 Important user information

Introduction

This section contains important user information about the product and this manual.

Read this before operating the product



All users must read the entire *Operating Instructions* before installing, operating, or maintaining the product.

Always keep the *Operating Instructions* at hand when operating the product.

Do not install, operate, or perform maintenance on the product in any other way than described in the user documentation. If you do, you may be exposed or expose others to hazards that can lead to personal injury and you may cause damage to the equipment.

Intended use of the product

The ÄKTA go instrument is intended for the purification of bio-molecules, in particular proteins, for research purposes. It is intended to be used by trained laboratory staff members in research laboratories within academia and industry.

The ÄKTA go instrument must not be used in any clinical procedures, or for diagnostic purposes.

System definition

In this manual, the combination of the ÄKTA go instrument and the UNICORN™ software is referred to as the system.

The ÄKTA go instrument without the software is referred to as the instrument.

Prerequisites

In order to operate ÄKTA go in the way it is intended:

- The user must know how to use a computer with the Windows operating system.
- The user must understand the concepts of liquid chromatography.
- The user must read and understand the *Safety instructions* chapter in the *Operating Instructions*.
- The ÄKTA go instrument must be installed in accordance with the site requirements and instructions in the *Operating Instructions*.

1.2 About this manual

Introduction

This section contains information about the purpose and scope of this manual, notes and tips, and typographical conventions.

Purpose of this manual

This manual provides information needed to install, operate and maintain the product in a safe way.

Scope of this manual

The *Operating Instructions* is valid for the ÄKTA go system.

Notes and tips

Note: *A note is used to indicate information that is important for trouble-free and optimal use of the product.*

Tip: *A tip contains useful information that can improve or optimize your procedures.*

Typographical conventions

Software items are identified in the text by ***bold italic*** text.

Hardware items are identified in the text by **bold** text.

Text that the user must either type exactly as shown in the manual, or that the software displays as a response (not a regular part of the graphic user interface), is shown by a monospaced typeface (for example, `Recipe Information`).

Tip: *The text can include clickable hyperlinks to reference information.*

1.3 Associated documentation

Introduction

This section describes the user documentation delivered with the product, and how to find related literature that can be downloaded or ordered from Cytiva.

ÄKTA go user documentation

The user documentation listed in the table below is available in printed format or as PDF file at [cytiva.com/aktago](https://www.cytiva.com/aktago) under **Related Documents**.

Translations of the original instructions are given in several languages and are contained in the CD provided with this manual, or can be found online at [cytiva.com/aktago](https://www.cytiva.com/aktago).

Documentation	Main contents
<i>ÄKTA go Unpacking Instructions (29383543)</i>	Information needed to handle the delivery package and unpack the ÄKTA go instrument.
<i>ÄKTA go Operating Instructions</i> (This instructions)	Information needed to install, operate, and maintain the ÄKTA go system in a safe way. Translations of the original instructions are given in several languages.
<i>ÄKTA go Cue Cards (29383545)</i>	Essential information to be kept near the ÄKTA go system.
<i>ÄKTA go User Manual (29391392)</i>	Additional detailed information on the system, component functions, and maintenance. Tips on how to get the most out of the system when running it.
<i>ÄKTA go Product Documentation (29434983)</i>	General specifications and list of materials in the flow path.
<i>ÄKTA avant, ÄKTA pure ÄKTA go, and ÄKTA pcc, Privacy and Security Manual, 29488174</i>	Describes the privacy and security considerations of the use of the system. The manual describes the expected intended use of the system, the privacy and security capabilities included, and how these capabilities are configured.
<i>ÄKTA avant, ÄKTA go, ÄKTA pcc, and ÄKTA pure Site Preparation Guide (29117084)</i>	Instructions on how to prepare the installation site for the ÄKTA go system.

UNICORN user documentation

The UNICORN user documentation is listed in the following table. It is available from the UNICORN software using the **Help** drop down menu **Contextual Help** under the heading **UNICORN contextual help and documentation**. It also can be download from cytiva.com/UNICORN under **Related Documents**.

Documentation	Main contents
<i>UNICORN Quick Installation Guide</i> ¹	Detailed instructions on how to install UNICORN.
<i>UNICORN Administration and Technical Manual</i> ²	<ul style="list-style-type: none"> • Overview and detailed description of network setup and complete software installation. • Administration of UNICORN and the UNICORN database.
<i>UNICORN Method Manual</i> ²	<ul style="list-style-type: none"> • Overview and detailed descriptions of the method creation features in UNICORN. • Workflow descriptions for common operations.
<i>UNICORN System Control Manual</i> ²	<ul style="list-style-type: none"> • Overview and detailed description of the system control features in UNICORN. • Includes general operation, system settings and instructions on how to perform a run.
<i>UNICORN Evaluation Manual</i> ²	<ul style="list-style-type: none"> • Overview and detailed descriptions of the Evaluation Classic³ module in UNICORN. • Description of the evaluation algorithms used in UNICORN.
Getting started with Evaluation (accessed through help in the UNICORN Evaluation module)	<ul style="list-style-type: none"> • Video clips showing common workflows in the Evaluation module. • Overview of features of the Evaluation module.
UNICORN Help	By pressing F1 , descriptions are displayed for the currently active pane or dialog box.

¹ The UNICORN Quick Installation Guide can be downloaded from cytiva.com/aktago.

² The current UNICORN version is added to the title of the manual.

³ Evaluation Classic is an advanced evaluation module that requires an extra license to run.

Additional literature

For practical tips on chromatography, refer to the handbooks available at cytiva.com/handbooks.

1.4 Abbreviations

Abbreviations

Abbreviation	Definition (English)	Translation (local language)
CIP	Cleaning In Place	Cleaning In Place
I/O box	Input/Output box	Input/Output box
RTU	Real-Time Unit	Real-Time Unit
UPS	Uninterruptible power supply	Uninterruptible power supply

2 Safety instructions

About this chapter

This chapter describes safety precautions, labels and symbols that are attached to the system. In addition, the chapter describes emergency and recovery procedures.

In this chapter

Section	See page
2.1 Safety precautions	11
2.2 Labels and symbols	15
2.3 Emergency procedures	16

Important



WARNING

All users must read and understand the entire contents of this general safety chapter, and the specific safety precautions information in each subsequent chapter of this manual to become aware of the hazards involved.

2.1 Safety precautions

Introduction

ÄKTA go is powered by mains voltage and handles materials that can be hazardous.

Before installing, operating, or maintaining the system, you must be aware of the hazards described in this manual.

Definitions

This user documentation contains safety notices (WARNING, CAUTION, and NOTICE) concerning the safe use of the product. See definitions below.



WARNING

WARNING indicates a hazardous situation which, if not avoided, could result in death or serious injury. It is important not to proceed until all stated conditions are met and clearly understood.



CAUTION

CAUTION indicates a hazardous situation which, if not avoided, could result in minor or moderate injury. It is important not to proceed until all stated conditions are met and clearly understood.



NOTICE

NOTICE indicates instructions that must be followed to avoid damage to the product or other equipment.

General precautions

The following general precautions must be considered at all times. There are also context related precautions, which are written in their respective chapters.



WARNING

Do not operate the product in any other way than described in the ÄKTA go user documentation.

**WARNING**

Only properly trained personnel are allowed to operate and maintain the product.

**WARNING**

Do not use any accessories not supplied or recommended by Cytiva.

**WARNING**

Do not use ÄKTA go if it is not working properly, or if it has suffered any damage, for example:

- damage to the power cord or its plug
- damage caused by dropping the equipment
- damage caused by splashing liquid onto it

**WARNING**

Access to power plug. Do not block access to the power outlet and power plug. The power cord with plug must always be easy to disconnect.

**WARNING**

In the event of a large spillage, disconnect the power cord from the wall socket.

**WARNING**

Always use appropriate Personal Protective Equipment (PPE) during operation and maintenance of this product.

**WARNING**

Hazardous substances and biological agents. When using hazardous chemical and biological agents, take all suitable protective measures, such as wearing protective clothing, glasses and gloves resistant to the substances used. Follow local and/or national regulations for safe operation and maintenance of this product.

**WARNING**

Spread of biological agents. The operator must take all necessary actions to avoid spreading hazardous biological agents. The facility must comply with the national code of practice for biosafety.

**WARNING**

High pressure. The product operates under high pressure. Wear protective glasses and other required Personal Protective Equipment (PPE) at all times.

Flammable liquids

**WARNING**

Explosion hazard: The product is **not approved** for use in a potentially explosive atmosphere. The product does not fulfill the requirements of the ATEX Directive.

**WARNING**

Explosion hazard. To avoid building up an explosive atmosphere when using flammable liquids, make sure that the room ventilation meets the local requirements.



WARNING

Explosion hazard due to leakage of flammable liquid. Make sure there is no leakage in the flow path of the instrument during the cleaning with a flammable liquid, such as ethanol. Carefully inspect the flow path for leakage, including the waste tubing and tighten any connectors if necessary.

2.2 Labels and symbols

Introduction

This section describes the nameplate, labels, and other safety and regulatory information attached to the product.

Nameplate

The nameplate provides information about the model, manufacturer, and technical data.

Precautions




WARNING

Access to power switch and power cord. The power switch must always be easy to access. The power cord must always be easy to disconnect.

Description of symbols and text

The following safety symbols and information may be present on the system label.

Label	Meaning
	Warning! Read the user documentation before using the system. Do not open any covers or replace parts unless specifically stated in the user documentation.
Voltage	Electrical rating: Voltage (VAC ~)
Frequency	Electrical rating: Frequency (Hz)
Max. Power	Electrical rating: Maximum power consumption (VA)
Protection Class	Degree of protection provided by the enclosure
Mfg. Date	Year (YYYY) and month (MM) of manufacture

2.3 Emergency procedures

Introduction

This section describes how to shut down the ÄKTA go instrument in an emergency situation, and the procedure for restarting the system.

The section also describes the result in the event of power failure.

Emergency shutdown

To shut down the instrument in an emergency, disconnect the instrument power cord from its power source. The power source can be an ordinary power outlet, or an UPS (Uninterruptible Power Supply) unit.

Any ongoing activity will be terminated immediately. Run data up to the time of the interruption will be saved.



NOTICE

Do not leave the instrument in an emergency stop condition. Flush the flow path with water or buffer when the emergency has been resolved.

Power failure

The result of a power failure depends on whether the system is equipped with a Real-Time Unit and whether the instrument or the computer is affected.

If power is lost to the ÄKTA go instrument, with or without a Real-Time Unit, the run is interrupted immediately. Run data collected up to the time of the power failure is saved in the UNICORN software.

If power is lost to the computer and the system is not equipped with a Real-Time Unit, the run is interrupted immediately. Run data collected up to the time of the power failure is saved in the UNICORN software.

If power is lost to the computer and the system is equipped with a Real-Time Unit, the run continues to completion and the run data is stored in the Real-Time Unit and uploaded to the computer once it is reconnected.

Note: *Connecting the instrument and computer to an uninterruptible power supply (UPS) can help to prevent loss of data and material during a power failure.*

Restart after emergency shutdown or power failure

Follow the steps below to restart the system after an emergency shutdown or power failure.

Step	Action
1	Reconnect the power cord.
2	Start the instrument by pressing the On/Off button on the instrument control panel.
3	Start the computer and the UNICORN software.
4	Re-establish connection between UNICORN and the instrument.
5	If the run has been aborted, recover or discard remaining sample and flush the flow path as appropriate.

3 System description

About this chapter

This chapter gives an overview of the ÄKTA go instrument and the UNICORN software.

In this chapter

Section		See page
3.1	ÄKTA go	19
3.2	Available modules	25
3.3	UNICORN	28

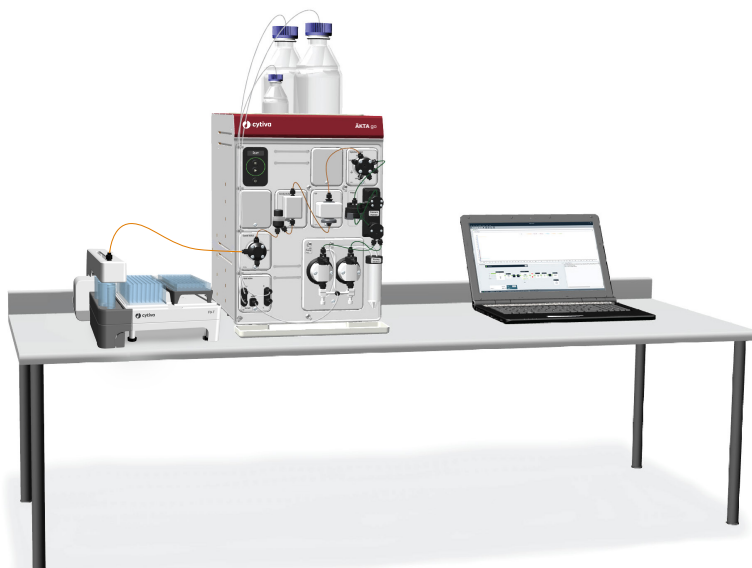
3.1 ÄKTA go

Introduction

This section provides an overview of the ÄKTA go instrument.

Illustration of the ÄKTA go system

The illustration below shows the ÄKTA go system. The computer with the UNICORN software is located on the right hand side of the instrument to make room for accessories on the left hand side of the instrument.

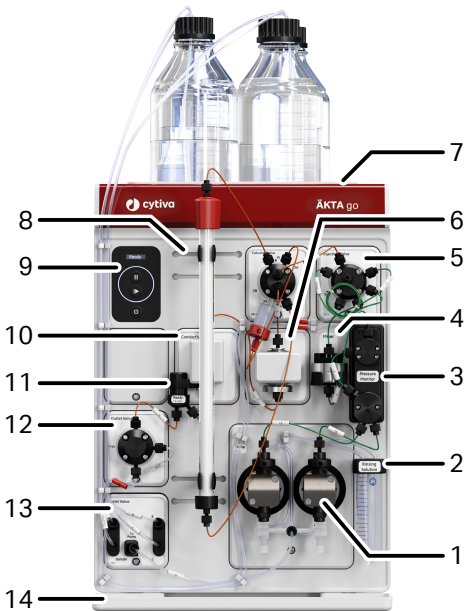


Exterior design

The liquid handling modules are placed on the front of the instrument. The instrument is equipped with trays to collect spillage, and adjustable feet to level the instrument. Buffer vessels can be placed on top of the instrument. Air ventilation, power cables, and data cables are located at the rear of the instrument.

Illustration of the ÄKTA go instrument

The illustration below shows the ÄKTA go instrument with all standard components denoted. The instrument in the illustration also has an optional column valve and a column connected.

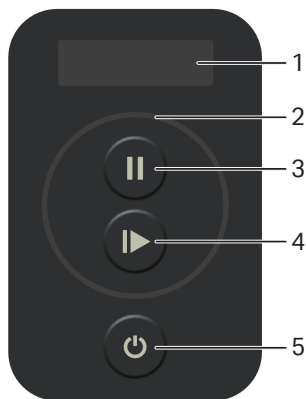


Part	Function
1	Pump
2	Pump rinsing solution tube
3	Pressure monitor
4	Mixer
5	Injection valve
6	UV monitor
7	Top tray
8	Holder rails
9	Instrument control panel
10	Conductivity monitor
11	Flow restrictor
12	Outlet valve
13	Inlet valve
14	Bottom tray

Illustration of the instrument control panel

The instrument control panel is located to the upper left on the front of the instrument. It shows current system status using LED light and status text. The pause and continue buttons can be used to control an ongoing run.

Note: Control panel buttons can be locked using the UNICORN software.





Part	Function
1	Display
2	Status indicator
3	Pause button
4	Run/Continue button
5	On/Off button

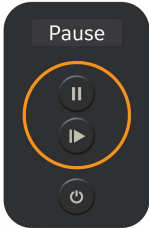
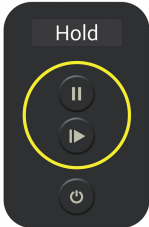
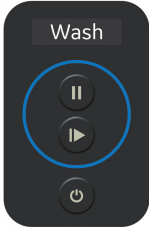
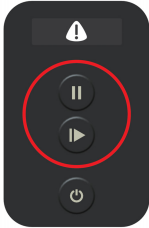
Status indicators



The display and status indicators on the instrument control panel indicate the current status of ÄKTA go.

The table below describes the different states that can be displayed.

State	Display	Description
Off		The instrument is turned off.
Turning on/off		The On/Off button is pressed and the instrument is turning on or off.

State	Display	Description
Offline	<p>The display toggles between Offline, the instrument IP address, and the Instrument serial number.</p>	Power is on, but the instrument has no communication with the UNICORN instrument server.
Connecting	<p>The status indicator flashes a white light.</p>	The instrument is connecting to the UNICORN instrument server.
Ready	<p></p>	The instrument is ready to use.
Run	<p></p>	A run is ongoing.

State	Display	Description
Pause		A run has been paused (pump is stopped).
Hold		A run has been put on hold (pump is still pumping at an unchanged flow rate).
Wash		A system wash or a pump wash instruction is ongoing.
Alarms and errors	 <p>The status indicator flashes a red light.</p>	The instrument has been paused due to an alarm or an instrument error. After investigating the cause of the error, acknowledge the alarm and continue the run in UNICORN.

State	Display	Description
Power-save	 <p>The status indicator is half-lit with a white light, flashing slowly.</p>	The instrument is in power-save mode.
Re-programming	 <p>Program</p>	A module is being re-programmed during an instrument configuration installation.

3.2 Available modules

Introduction

The ÄKTA go instrument is delivered with standard modules installed. There is space for installation of six additional modules in the system, two inside the chassis and four connected via cables at the rear of the instrument. This section describes the standard and optional modules.

Standard modules

The following modules are delivered installed in the instrument.

Module	Description
Inlet valve K9	An inlet valve for buffers, sample, and cleaning solution. Creates a gradient by switching between inlets A and B .
Pump P9-S	A high precision pump that delivers buffer or sample.
Pressure monitor R9	A pressure monitor that measures the pressure directly after the pump.
Mixer	A 1 mL static mixer that mixes the buffers delivered by the pump.
Injection valve V9-J	An injection valve that injects sample onto the column.
UV monitor U9-L	A LED UV monitor that measures the UV absorbance at 280 nm of buffers and eluted proteins. Includes a UV cell with a path length of 2 mm. An optional UV cell with a path length of 5 mm is available. The LED lamp does not require warming before use and does not heat the sample.
Conductivity monitor C9	A conductivity monitor that measures the conductivity of buffers and eluted proteins.
Outlet valve V9-Os	An outlet valve that directs the flow from the instrument to a fraction collector, outlet 1, or to waste.

Optional modules

The following modules can be added to the system. A maximum of six optional modules can be connected to the system.

Module	Description
External air sensor L9-1.5	An air sensor used to either complete sample loading via the pump, or to detect if the system has run out of buffer, depending on placement. Tubing connector: 5/16" + Ferrule (yellow), 1/8".
External air sensor L9-1.2	An air sensor used to complete sample loading via the pump. Do not place L9-1.2 between the inlet valve K9 and the pump. Tubing connector: Union 1/16" Male, 5/16" Female fitting tubing connector, 5/16" + Ferrule (yellow), 1/8". Do not connect this air sensor with narrow tubing (o.d. 1/16") because the high flow rate during a pump wash might cause cavitation.
Inlet valve V9-lmA	An inlet valve that enables six buffer inlets.
Inlet valve V9-lmB	An inlet valve that enables six buffer inlets.
Inlet valve V9-lmS	An inlet valve that enables five sample inlets and one buffer inlet.
Column valve V9-Cm	A column valve that can connect up to three columns to the instrument. Directs the flow onto one column at a time and allows for flow in two directions (upflow and downflow).
Column valve V9-C	An advanced column valve that can connect up to five columns to the instrument. Directs the flow onto one column at a time and allows for flow in two directions (upflow and downflow). This module contains two pressure sensors that enable pre-column pressure and delta-column pressure signals.
pH valve V9-pH	A pH valve that enables in-line monitoring of pH during the run.
Outlet valve V9-O	An outlet valve that directs the flow from the instrument to a fraction collector, to any of the ten outlet ports, or to waste.
Fraction collector F9-R	A fraction collector with up to 175 fractions.
Fraction collector F9-T	A fraction collector to fractionate in two plates, or small tubes placed in racks, and in four 50 mL tubes.
I/O-box E9	An I/O box that sends and receives analog or digital signals to and from external equipment.

Note: *You must create new methods if you add or remove modules from the system, since the available software instructions are updated when modules are added or removed.*

Real-Time Unit

The system can be equipped with a Real-Time Unit (RTU), which must be installed by a Cytiva service representative. The RTU can be used in certain network environments to make sure the run continues if the computer is rebooted or otherwise locked due to, for example, software updates. The method is automatically downloaded to the RTU when the method is started. Instructions are sent from the RTU to the instrument and data are collected in the RTU during the run. The result is uploaded from the RTU to the computer when the connection is restored.



NOTICE

Damage to RTU. To shut down the system, make sure to press the On/Off button for a short time. When the button is pressed for an extended period of time, the system is promptly shut down, which may damage the Real-Time Unit (RTU).

3.3 UNICORN

Introduction

ÄKTA go systems are controlled by UNICORN software running on an external computer. This section gives a brief overview of the UNICORN modules and the **System Control** user interface. Refer to UNICORN user documentation for more information.

The examples given in this manual are from UNICORN 7.9.

UNICORN modules overview

UNICORN consists of four modules: **Administration**, **Method Editor**, **System Control**, and **Evaluation**. The main functions of each software module are described in the following table.

Software module	Main functions
Administration	Perform user and system setup, system log, and data-base administration.
Method Editor	Create and edit methods using one or a combination of: <ul style="list-style-type: none"> • Predefined methods with built-in application support • Drag-and-drop function to build methods with relevant steps • Method text editing
System Control	Start, monitor, and control runs. The current flow path is illustrated in the Process Picture , which allows manual interactions with the system and provides feedback on run parameters.
Evaluation	Open results and evaluate runs. The Evaluation module includes a user interface optimized for workflows such as quick evaluation, comparing results, and working with peaks and fractions. Advanced features requires Evaluation Classic , available from Cytiva.

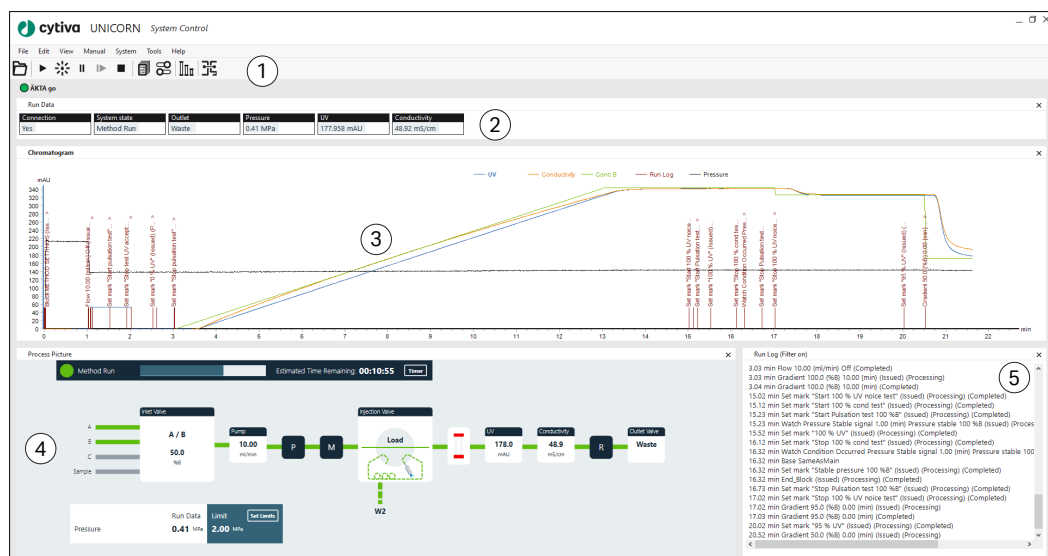
When working with the software modules **Administration**, **Method Editor**, **System Control**, and **Evaluation Classic** it is possible to access descriptions of the active window or software instruction by pressing the **F1** key. The module **Evaluation** contains a guide with video clips.

Opening a UNICORN module

Modules to open are selected at log in, but can also be opened from another module when the software is already open. In the **Administration**, **Method Editor**, or **System Control** modules, to open a software module, click **Tools** and select the applicable module. When in the **Evaluation** module, to open a software module, click **File** → **Applications** and select the applicable module.

Illustration of the System Control user interface











The illustration below shows the **System Control** module.



Part	Description
1	The toolbar buttons are used for quick access to instrument controls. For descriptions, see <i>System Control toolbar buttons</i> below.
2	The run data field shows the value of run data in boxes. This field is hidden by default. To make this field visible, go to View → run data .
3	The Chromatogram pane illustrates the chromatogram of the run.
4	The interactive Process Picture pane allows manual interactions with the system, illustrates the current flow path, and provides feedback on component status and run parameters.
5	The Run Log pane shows all registered events during the run.

System Control toolbar buttons

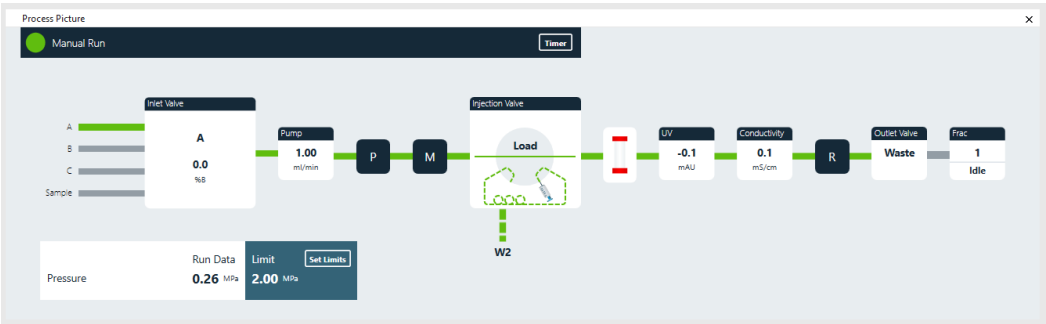
The following table shows the **System Control** toolbar buttons.

Button	Function	Button	Function
	Open Method Navigator This button opens the Method Navigator where saved methods are listed.		Run This button starts a method run. The last method run will be started.
	Hold This button suspends the method run, while current flow rate and valve positions are sustained.		Pause This button suspends the method run and stops all pumps.
	Continue This button resumes a held or paused method run.		End This button permanently ends the run.
	Documentation This button opens a dialog containing information about the system and the current run.		Customize This button opens the Customize dialog box where curve settings, run data groups and run log contents can be set.
	Column Handling This button opens the Column Handling tool, which contains a column list, with parameters for Cytiva columns. With an additional license, a Column Logbook to keep track of user-purchased columns is also available.		Connect to Systems This button opens the Connect to Systems dialog box where systems can be connected, and currently connected users are displayed.

Process Picture pane

The most commonly used manual interactions can be executed using the **Process Picture**. Click on the different parts of the **Process Picture** pane to interact with the system.

For a complete list of manual instructions, go to **Manual** → **Execute manual instructions**.



4 Installation

About this chapter

This chapter provides the instructions necessary to enable users to install the instrument and the software. Read the entire Installation chapter before starting to install ÄKTA go.

Note: For information how to unpack the ÄKTA go instrument and how to lift the instrument onto a laboratory bench refer to the ÄKTA go Unpacking instructions.

In this chapter

Section	See page
4.1 Site preparation	33
4.2 Hardware installation	44
4.3 Software installation	54
4.4 Start UNICORN and connect to system	59

4.1 Site preparation

Introduction

This section describes the site planning and the preparations necessary for the installation of ÄKTA go.

The performance specifications of the system can be met only if the laboratory environment fulfills the requirements stated in this chapter.

In this section

Section		See page
4.1.1	Delivery, storage, and unpacking	34
4.1.2	Space requirements	36
4.1.3	Site environment	39
4.1.4	Power requirements	41
4.1.5	Computer requirements	42
4.1.6	Required materials	43

4.1.1 Delivery, storage, and unpacking

ÄKTA go Unpacking Instructions



CAUTION

Heavy object. Use suitable precautions when moving the instrument. Two people are recommended to lift the instrument safely.

For information on how to receive and store the delivery box, and to unpack the ÄKTA go instrument refer to the *ÄKTA go Unpacking Instructions* attached on the delivery box. It is also available on the web. Two people are required to safely unpack the instrument, and no special equipment is needed.

When you receive the delivery

- Record on the receiving documents if there is any apparent damage on the delivery box. Inform your Cytiva representative of such damage.
- Move the delivery box to a protected location indoors.

Delivery box

The ÄKTA go instruments are shipped in a delivery box with the following dimensions and weight:

Contents	Dimensions (mm) w × h × d	Weight
ÄKTA go instrument with accessories	509 × 566 × 774	39 kg

Storage requirements

The delivery box should be stored in a protected place indoors. The following storage requirements must be fulfilled for the unopened box:

Parameter	Allowed range
Ambient temperature, storage	-25°C to 60°C for 48 h
Relative humidity	up to 90% atmospheric humidity at 40°C

Moving the ÄKTA go instrument

The instrument is heavier at the front. Do not tip the instrument when lifting. The illustration below shows the recommended way to lift the ÄKTA go instrument.



4.1.2 Space requirements

Introduction

This section describes the requirements for the laboratory bench on which the instrument is placed.

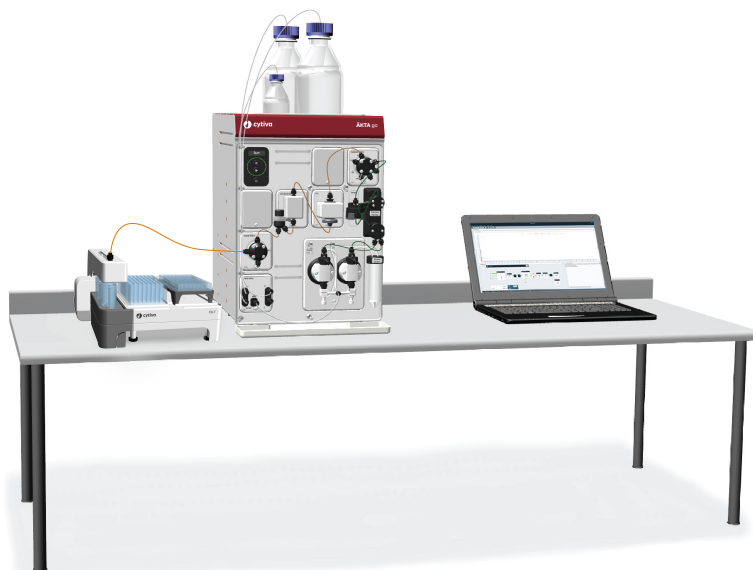


WARNING

Access to power plug. Do not block access to the power outlet and power plug. The power cord with plug must always be easy to disconnect.

Laboratory bench

The bench must be clean, flat and stable, and of sufficient size for the ÄKTA go system with accessories. The computer is located on the right side of the instrument to make room for accessories on the left side of the instrument.



See the table below for a summary of space requirement.

Parameters	Specifications
Min. length of the laboratory bench (with one fraction collector, without the computer)	76 cm
Min. depth of the laboratory bench	60 cm

Parameters	Specifications
Min. height above the laboratory bench	48 cm
Space for handling of buffers and samples	10 cm
Free space around the instrument to allow for adequate ventilation	10 cm
Location of the computer with respect to instrument	Right

Dimensions and weight



Parameter	Value
W (width)	335 mm
H (height)	482 mm
D (depth)	464 mm

Parameter	Value
Weight	27 kg
Allowed weight of buffer bottles on the top tray	10 kg

4.1.3 Site environment

Introduction

This section describes the environmental requirements for installation of ÄKTA go.

Environmental requirements

The following general requirements must be fulfilled:

- The room must have exhaust ventilation.
- The instrument must not be exposed to direct sunlight.
- Dust in the atmosphere must be kept to a minimum.

The installation site must comply with the following specifications.

Parameter	Requirement
Allowed location	Indoor use only
Ambient temperature, operating	4°C to 35°C
Ambient temperature, storage	-25°C to 60°C for 48 h
Relative humidity	20% to 95%, non-condensing
Altitude, operating	Up to 2000 m ¹
Pollution degree of the intended environment	Pollution degree 2

¹ If the installation is situated at a higher altitude, the customer is responsible for establishing that the ÄKTA go instrument can be used safely in accordance with local regulations.

Instrument ventilation

There should be at least 10 cm clearance at the back of the instrument to allow adequate air circulation.



NOTICE

Do not block the air vents on the rear of the instrument.

Heat output

The heat output is listed in the table below.

Component	Heat output
ÄKTA go instrument	Typical 100 W Max 150 W Power save < 20 W

For heat output of the computer, refer to the manufacturer's specifications.

4.1.4 Power requirements

Introduction

This section describes the power supply requirements for ÄKTA go.

Electrical power requirements



WARNING

Protective ground. The product must always be connected to a grounded power outlet.



WARNING

Power cord. Only use power cords with approved plugs delivered or approved by Cytiva.

The following table specifies the power requirements for the ÄKTA go instrument. For power requirements for the computer, refer to the manufacturer's specifications.

Parameter	Requirement
Supply voltage	100 to 240 V~
Frequency	50/60 Hz
Transient overvoltages	Overvoltage category II
Max power consumption	300 VA

4.1.5 Computer requirements

Introduction

This section describes the computer requirements for the instrument.

General requirements

ÄKTA go instruments are controlled by UNICORN software running on an external computer. The computer is not included with the ÄKTA go instrument.

A suitable computer may be ordered from Cytiva or obtained from a third party supplier.



WARNING

Fire hazard. Make sure the installed computer and other equipment has appropriate protection for the intended environment. Exposure to liquids or moisture can cause a short circuit and fire.



NOTICE

Any computer used with the equipment must comply with IEC 60950 and be installed and used according to the manufacturer's instructions.

For information on computer specifications, see [cytiva.com/UNICORNPCspecifications](https://www.cytiva.com/UNICORNPCspecifications).

For information about compatibility between UNICORN versions, the supported operating systems, database versions, and instrument configuration, see the UNICORN compatibility matrix at [cytiva.com/UNICORNcompatibility](https://www.cytiva.com/UNICORNcompatibility).

Network connection requirements

If the computer is to be connected to a network by a network cable, two Ethernet ports are required on the computer, one for connection to the ÄKTA go instrument and one for connection to a network.

Network settings are described in UNICORN user documentation.

4.1.6 Required materials

Introduction

This section describes the materials required for the installation and operation of the ÄKTA go instrument.

Solutions

The solutions listed in the following table are required during the installation procedure and should be provided at the installation site.

Buffer/solution	Required volume	Scope of use
Distilled water	1000 mL	System test, Pump test, and Mixer test If applicable, Fraction Collector F9-R test, Fraction Collector F9-T test and/or Column Valve V9-C test
1.0% acetone and 1.0 M NaCl in distilled water	1000 mL	System test and Mixer test
20% ethanol in distilled water	1000 mL	Priming of the pump rinsing system

Note: *The solutions must be freshly prepared, in order for the installation procedure to be successful. Delivery and installation might not take place at same time.*

Note: *Make sure that the solutions are at the same temperature as the intended location of the ÄKTA instrument.*

4.2 Hardware installation

Introduction

This section describes the installation procedures for ÄKTA go.

In this section

Section		See page
4.2.1	Connect the system	45
4.2.2	Install waste tubing	48
4.2.3	Prepare the pump rinsing system	50
4.2.4	Start the instrument	53

4.2.1 Connect the system

Introduction

The following connections must be made:

- Power supply to the ÄKTA go instrument and the computer
- Network connection between the computer and the ÄKTA go instrument



WARNING

Access to power plug. Do not block access to the power outlet and power plug. The power cord with plug must always be easy to disconnect.



WARNING

Protective ground. The product must always be connected to a grounded power outlet.



WARNING

Power cord. Only use power cords with approved plugs delivered or approved by Cytiva.

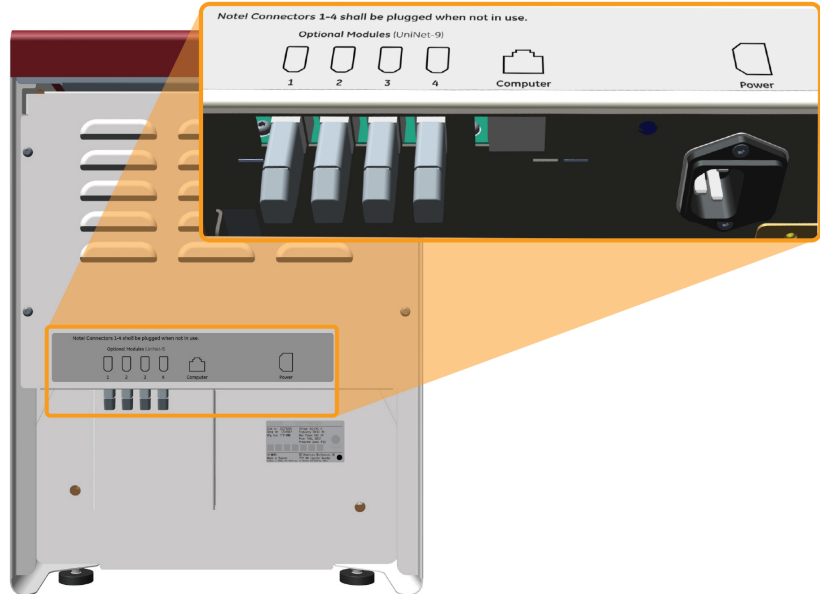


WARNING

Supply voltage. Before connecting the power cord, make sure that the supply voltage at the wall outlet corresponds to the marking on the instrument.

Connector illustration

The illustration below shows where the connectors are located on the ÄKTA go instrument. For connectors on the computer equipment, refer to the manufacturer's documentation.




Connect to power

Follow the steps below to connect power to the ÄKTA go instrument and the computer.

Step	Action
1	<p>Select the correct power cord to be used. Each instrument is delivered with 2 alternative power cords:</p> <ul style="list-style-type: none"> • Power cord with US-plug, 2 m • Power cord with EU-plug, 2 m <p>Recycle the power cord that is not to be used.</p>
2	<p>Connect the power cord to the Power input connector on the back of the instrument and to a grounded power outlet 100 to 240 V~, 50/60 Hz.</p>
3	<p>Connect the computer to a power source using the manufacturer's instructions.</p>

Connect ÄKTA go to the computer

Follow the steps below to connect the ÄKTA go to the computer.

Step	Action
1	<p>Connect a network cable to the back of the instrument. The appropriate port at the back of the instrument is indicated by this symbol:</p>  <p>Computer</p>
2	<p>Connect the other end of the network cable to the appropriate connector on the computer.</p>
3	<p>Make sure that the IP address of the instrument is on the same subnet as the IP address of the port used in the computer.</p> <p>Note:</p> <p><i>The IP address of the instrument is displayed on the instrument control panel when the instrument is powered on but not connected to UNICORN.</i></p>
4	<p>If the computer is to be connected to a network, connect a network cable between the computer and a network wall outlet.</p>

4.2.2 Install waste tubing

Introduction

The table below lists the waste tubing of the instrument and where it is located. Make sure that the waste tubing is connected to the correct positions on the modules.

Module	Tubing connections	Location of tubing
Injection valve	Waste ports W1 and W2	Front of the ÄKTA go instrument.
Outlet valve (optional)	Waste port W	Front of the ÄKTA go instrument.
pH valve (optional)	Waste port W3	Front of the ÄKTA go instrument.

Prepare waste tubing

Follow the instructions below to prepare the waste tubing.



CAUTION

Fasten the waste tubing. During operation at high pressure the ÄKTA go instrument may release bursts of liquid in the waste tubing. Make sure that all waste tubing is securely fastened to the ÄKTA go instrument and to the waste vessel.



CAUTION

Make sure that the waste vessel will hold all the produced volume of the run. For ÄKTA go, a suitable waste vessel should typically have a volume of 2 to 10 liters.



CAUTION

Cut injuries. The tubing cutter is very sharp and must be handled with care to avoid injuries.



NOTICE

The highest level of liquid in the waste vessel must be lower than the Waste port **W** on the Outlet valve of the ÄKTA go instrument.

Step	Action
1	Insert the waste tubing from all installed modules in a suitable vessel.
2	Cut the waste tubing to appropriate length. It is important that the tubing is not bent.



Note:

If the tubing is too short, replace it with new tubing. Do not lengthen the tubing as this might cause obstruction of the tubing.

3	Fasten all waste tubing to the waste vessel.
---	--

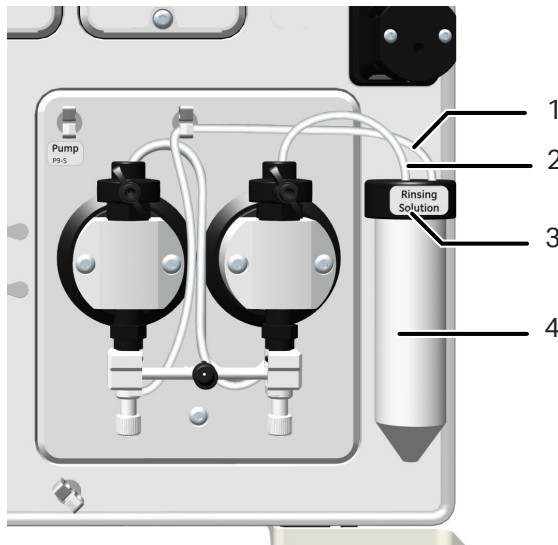
4.2.3 Prepare the pump rinsing system

Introduction

The pump rinsing system protects the pump seals from damage caused by precipitated buffer remaining in the system. The seal prevents leakage between the pump chamber and the drive mechanism of the pump.

Illustration of the pump rinsing system

The illustration below shows the parts and tubing of the pump rinsing system.



Part	Description
1	Inlet tubing
2	Outlet tubing
3	Rinsing solution tube holder
4	Rinsing solution tube

Prime the pump rinsing system

Follow the instructions below to fill the pump rinsing system with rinsing solution.

Step	Action
------	--------

- | | |
|---|---|
| 1 | Unscrew and remove the rinsing solution tube from the holder. |
|---|---|

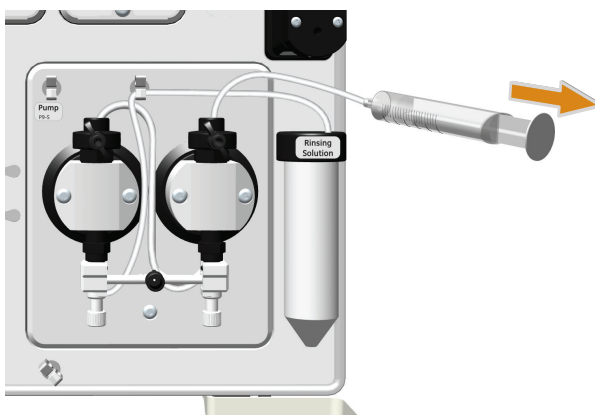


- | | |
|---|---|
| 2 | Fill the rinsing solution tube with 50 mL of 20% ethanol or aqueous buffer. |
| 3 | Screw the rinsing solution tube back into the holder. |
| 4 | Insert the inlet tubing into the solution in the rinsing solution tube. |

Note:

Make sure that the inlet tubing reaches close to the bottom of the rinsing liquid tube.

- | | |
|---|--|
| 5 | Connect a 25 to 30 mL syringe to the outlet tubing. Draw liquid slowly into the syringe until the rinsing system tubing is filled. |
|---|--|



- | | |
|---|--|
| 6 | Disconnect the syringe and discard its contents. |
| 7 | Insert the outlet tubing into the liquid in the rinsing solution tube. |

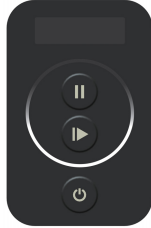
Step	Action
8	Unscrew the rinsing solution tube and fill it with 50 mL of 20% ethanol or aqueous buffer.

4.2.4 Start the instrument

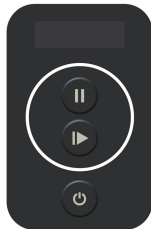
Follow the steps below to start the instrument.

Step	Action
------	--------

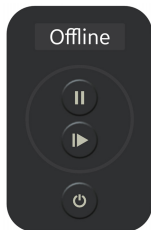
- | | |
|---|---|
| 1 | Turn on the instrument by pressing the On/Off button. |
|---|---|



- | | |
|---|--|
| 2 | The instrument control panel displays a white light for approximately two seconds. |
|---|--|



- | | |
|---|--|
| 3 | The instrument is on, but is not connected to the UNICORN instrument server. |
|---|--|



Result:

The displays toggles between **Offline**, the instrument IP address, and the instrument serial number.

4.3 Software installation

Introduction

This section provides an overview of how to install UNICORN and adapt the software to your instrument. For more information, refer to the *UNICORN Quick Installation Guide*.

The software should be installed by someone assigned to be a UNICORN system administrator at the site. Detailed information about software installation and configuration is available in the *UNICORN Administration and Technical Manual*.

In this section

Section		See page
4.3.1	Download and Install UNICORN	55
4.3.2	Download the Instrument Configuration	56
4.3.3	Adapt UNICORN to your system	57

4.3.1 Download and Install UNICORN

UNICORN is delivered via e-Delivery. An URL to the e-Delivery and Activation ID are delivered upon ordering the ÄKTA go system.

Follow the steps below to install the UNICORN software. For more information on installing UNICORN, Windows settings, and configuring the e-license, refer to the *UNICORN Quick Installation Guide*.

Step	Action
1	Download UNICORN from the e-Delivery portal.
2	Start the installation wizard.
3	On the welcome dialog box, click Next .
4	Select installation type, Full installation or Custom installation , and click Next . Note: <i>Any missing prerequisites will be installed. The computer might need to be restarted several times during the installation.</i>
5	Download an e-License from the e-Delivery portal and configure the e-license for the UNICORN installation.
6	Adapt UNICORN to your system, see Section 4.3.3 Adapt UNICORN to your system, on page 57 . Note: <i>The computer and the instrument must have IP addresses on the same subnet.</i>

4.3.2 Download the Instrument Configuration

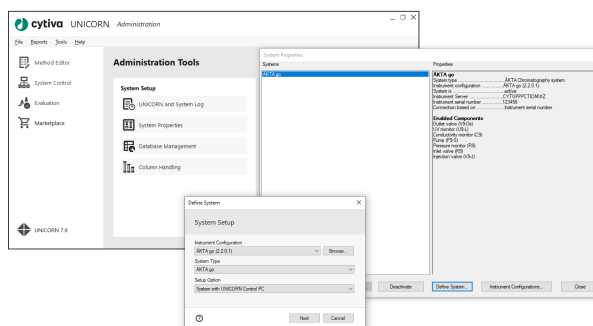
An instrument configuration is used to adapt UNICORN to your instrument. Follow the instructions below to import the Instrument Configuration into the UNICORN software.

Step	Action
1	Go to cytiva.com .
2	Click RELATED DOCUMENTS .
3	Click SOFTWARE .
4	Download the Instrument configuration software .
5	Use the downloaded instrument configuration to define your system, see Section 4.3.3 Adapt UNICORN to your system, on page 57 .

4.3.3 Adapt UNICORN to your system

To be able to connect to the system, UNICORN must be adapted to the instrument with the correct instrument configuration installed and the correct modules selected in the software. This is done following the steps below.

Step	Action
1	Download the latest instrument configuration for ÄKTA go from cytiva.com/aktago , see Section 4.3.2 Download the Instrument Configuration, on page 56 .
	Note: <i>An instrument configuration is used to adapt UNICORN to your instrument.</i>
2	In the Administration module, click System Properties → Instrument Configurations and import the downloaded instrument configuration.
3	Click Define System to define a new system with the imported instrument configuration.



Note: *A detailed description of how to adapt UNICORN to your system, including system setup, is found in the UNICORN Administration and Technical Manual.*

4.4 Start UNICORN and connect to system

Introduction

This section describes how to start and log on to UNICORN, and how to connect to the system in UNICORN.

Prerequisites


For UNICORN to be correctly installed, the following conditions must be set:


- the IP address of the computer must be set to the same subnet as that of the instrument
- the e-license needs to be downloaded and configured for the computer
- the system has to be defined with the correct modules in UNICORN

Start UNICORN and log on

Follow the instructions to start UNICORN and log on to the program.

Step	Action
1	Double-click the UNICORN icon on the desktop.
2	In the Log On dialog box, select User Name and enter Password .


UNICORN™ 7.9
Version 7.9.0.2440

☐ Use Windows Authentication 

User Name

Default

Password

Domain

Access

Administrators

☒ Administration

☒ System Control

☐ Method Editor

☒ Evaluation

LOG IN

CANCEL

3	Tick the boxes for the UNICORN modules that you want to start.
---	--

Step	Action
4	Click OK .

Connect to system

Follow the instructions to connect to the system in UNICORN.

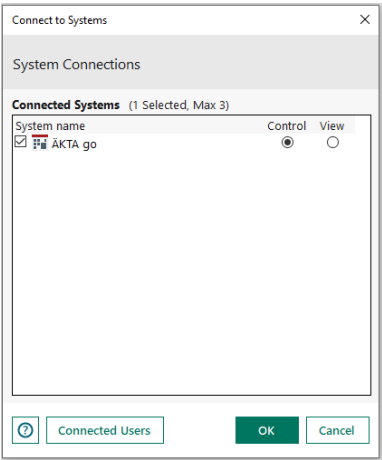
Note: *The system must have been defined by the UNICORN system administrator for it to be present in the database.*

Step	Action
1	In the System Control module, click the Connect to Systems button.



2 In the **Connect to Systems** dialog box:

- Select a system checkbox.
- Click **Control** for that system.
- Click **OK**.



Tip:
If UNICORN is unable to connect to the selected instrument, try restarting the computer. See [Chapter 8 Troubleshooting, on page 102](#) for more information.

Set up **Power-save**

To minimize power consumption when the system is not used, there is a **Power-save** function in UNICORN. Follow the steps below to activate the **Power-save** function.

Step	Action
------	--------

- | | |
|---|--|
| 1 | In System Control click System → Settings → Advanced . |
| 2 | Turn on the Power-save function. |
| 3 | Enter a Time for the Power-save |
| 4 | Click OK . |

5 Prepare the system for a run

About this chapter

This chapter gives instructions on how to prepare the ÄKTA go system for a run and what to do before the first run.

In this chapter

Section	See page
5.1 Prepare the flow path	64
5.2 Prime inlets and purge pump heads	66
5.3 Performance tests	72
5.4 Connect a column	75
5.5 Pressure alarms	78
5.6 Prepare for a run at low temperature	79

Safety Precautions



WARNING

Explosion hazard. To avoid building up an explosive atmosphere when using flammable liquids, make sure that the room ventilation meets the local requirements.



CAUTION

Fire hazard. Before starting the system, make sure that there is no leakage of flammable or potentially explosive liquids.



WARNING

Cutting risk. Use the protective jacket and do not exceed the pressure specifications of the Superloop. Overpressure can cause the glass in the Superloop to break.

**CAUTION**

Fasten bottles. Always fasten bottles to the rails at the front panel. Use appropriate holders for bottles. Shattered glass from falling bottles may cause injury. Spilled liquid may cause fire hazard and personal injury.

**CAUTION**

Max. weight on top tray. Do not place containers with a volume of more than 2 liters each on the top tray. The total allowed weight on the top tray is 10 kg.

**CAUTION**

Avoid spillage and overflow during collection. Make sure that:

- An outlet tubing is connected to each outlet port.
- A vessel is connected to collect from each outlet tubing.
- The fraction collector is connected and loaded with correct plates or tubes.

**CAUTION**

Avoid spillage and overflow from waste. Make sure that the waste tubing is inserted in an appropriate waste container and secured in place.

**CAUTION**

Cut injuries. The tubing cutter is very sharp and must be handled with care to avoid injuries.

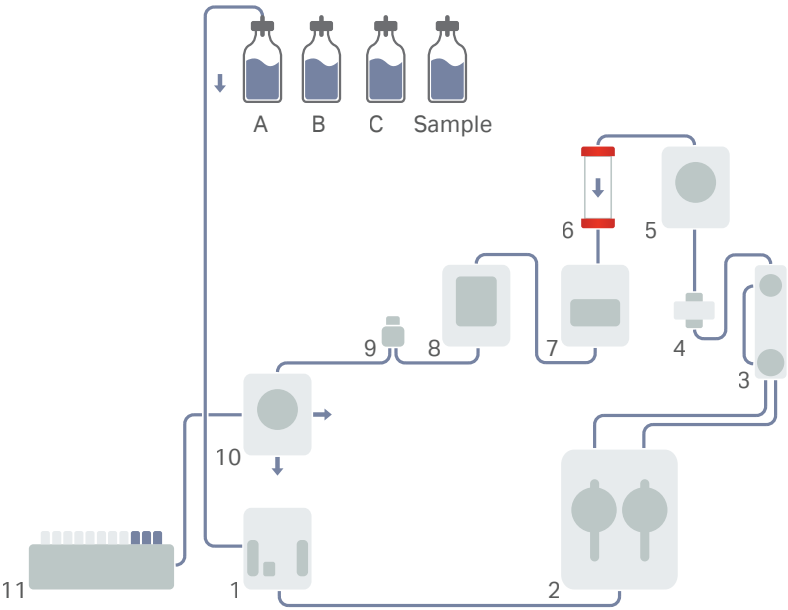
5.1 Prepare the flow path

Introduction

The ÄKTA go instrument, as delivered, is prepared with a default flow path. The modules in this flow path must be defined in the software, see [Section 4.3.3 Adapt UNICORN to your system, on page 57](#). It is possible to remove modules from the flow path and to add some additional valves and monitors. If the modules in the flow path on the instrument are changed, the **System Properties** in the software must be updated.

Illustration of the flow path

The illustration below shows the flow path for a standard configured ÄKTA go instrument with an optional fraction collector connected. The individual instrument modules are presented in the table below. The system configuration is defined by the user in the UNICORN **Administration** module.



Part	Description	Part	Description
1	Inlet valve K9	7	UV monitor U9-L
2	Pump P9-S	8	Conductivity monitor C9
3	Pressure monitor R9	9	Flow restrictor FR-902

Part	Description	Part	Description
4	Mixer	10	Outlet valve V9-Os/V9-O
5	Injection valve V9-J	11	Fraction collector F9-T/F9-R
6	Column		

Prepare the waste tubing

Make sure that the waste tubing is prepared according to the instructions in [Section 4.2.2 Install waste tubing, on page 48](#).

Prepare the outlet tubing

Connect tubing to the outlet ports of the outlet valve that are to be used during the run.

If no fraction collector is used, immerse the outlet tubing in a suitable flask.

If a fraction collector is used, make sure that 400 mm tubing (or 800 mm tubing if a tunnel is used) is connected between the fraction collector and the **Frac** port on the outlet valve, and prepare the fraction collector for a run. For more information, refer to the *Operating Instructions* of the fraction collector.

Plug unused valve ports

It is recommended to plug all unused valve ports with stop plugs before starting a run. Refer to the *ÄKTA go User Manual* for information about connectors.



CAUTION

Make sure there is a sample loop or stop plugs in place in the loop positions in the injection valve to avoid leakage during valve turns.

5.2 Prime inlets and purge pump heads

Introduction

Before using the pump, it is important to prime all inlets and purge the pump heads, that is, to fill the inlets and pump heads with liquid so that no air remains inside.

Overview

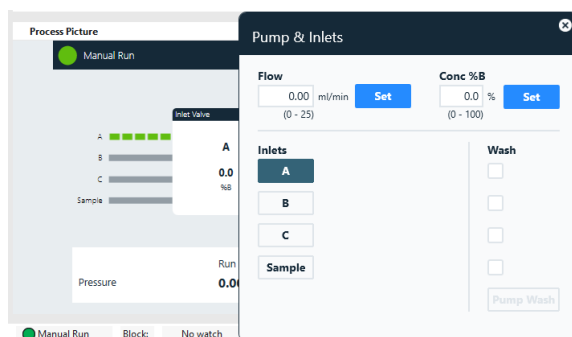
The procedure consists of the following stages:

1. Prime the inlet tubing
2. Purge the pump and confirm that it is free from air

Prime the inlet tubing

Follow the steps below to prime all inlet tubing that is to be used during the run.

Step	Action
1	Make sure that all inlet tubing that is to be used during the run is placed in the correct buffer.
2	Open the System Control module.
3	In the Process Picture pane, click on Inlet Valve and select one of the inlets to be primed.

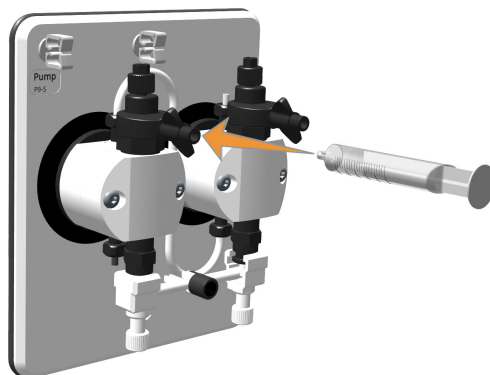


Result:

The inlet valve opens the selected inlet.

Step	Action
------	--------

- | | |
|---|--|
| 4 | Connect a 25 to 30 mL syringe to the purge valve of one of the pump heads. |
|---|--|



- | | |
|---|--|
| 5 | Open the purge valve by turning it counter-clockwise one and a half turns. Draw liquid slowly into the syringe until liquid reaches the pump and the inlet tubing is filled with liquid. |
| 6 | Close the purge valve by turning it clockwise. Disconnect the syringe and discard its contents. |
| 7 | Repeat steps 3 to 6 for each inlet that is to be used during the run. |

Purge the pump and confirm it is free from air

Follow the steps below to purge the two pump heads of the pump so that they are free from air.

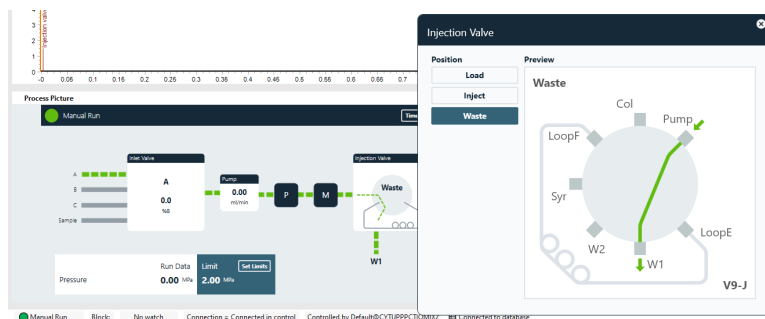
Step	Action
------	--------

- | | |
|---|--|
| 1 | In the Process Picture pane, click Injection valve and select Waste . |
|---|--|

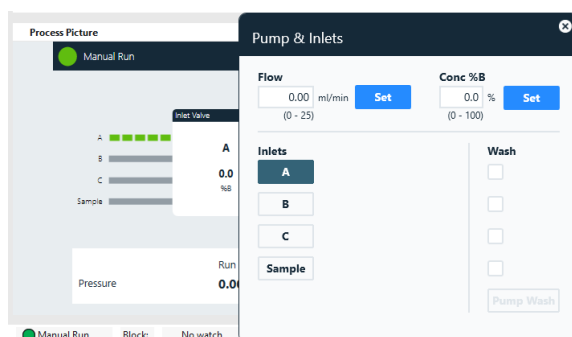
Step	Action
------	--------

Result:

The injection valve switches to waste position and opens inlet **A**.



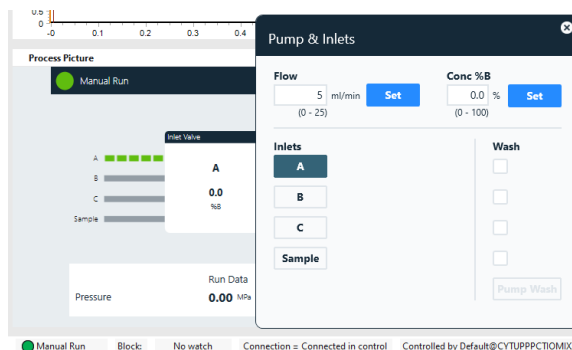
- 2 In the **Process Picture** pane, click **Inlet Valve** and select the inlet that will be used at the beginning of the run.



Result:

The inlet valve opens the selected inlet.

- 3 Under **Flow**, enter 5 mL/min and click **Set**.

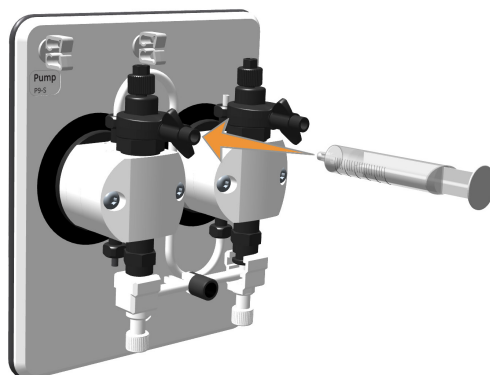


Step	Action
------	--------

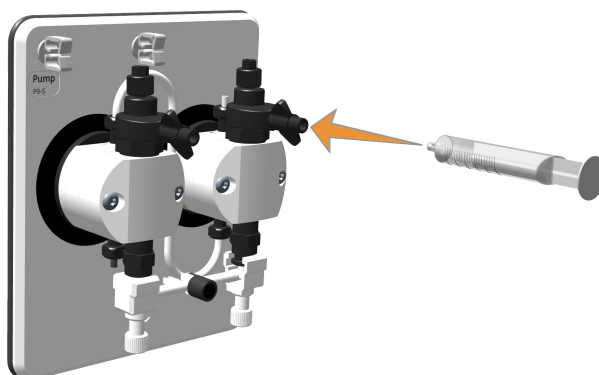
Result:

A system flow starts.

- | | |
|---|---|
| 4 | Connect a 25 to 30 mL syringe to the purge valve of the left pump head. |
|---|---|



- | | |
|---|--|
| 5 | Open the purge valve by turning it counter-clockwise about one and a half turns. Draw liquid slowly into the syringe, at a rate of approximately 1 mL per second, until there are no air bubbles in the liquid that reaches the syringe. |
| 6 | Close the purge valve by turning it clockwise. Disconnect the syringe and discard its contents. |
| 7 | Keep the system flow running. Connect the syringe to the purge valve of the right pump head, and repeat steps 5 and 6. |



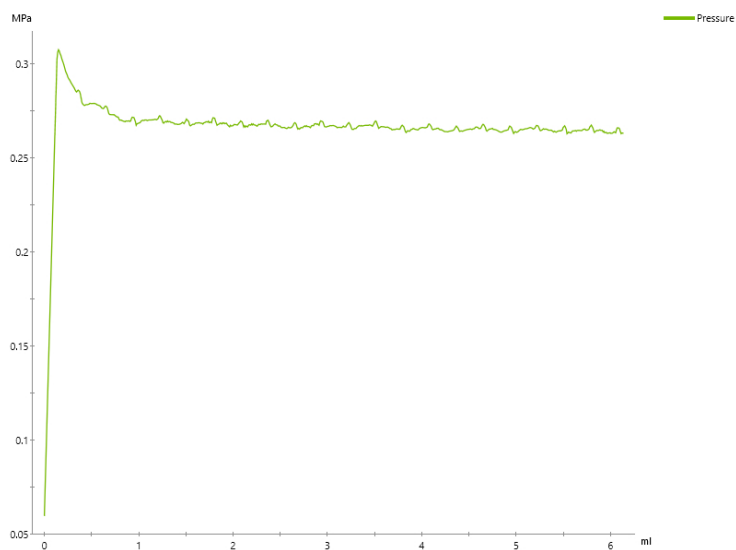
Step	Action
------	--------

- | | |
|---|--|
| 8 | In the Chromatogram pane, check the pressure curve. If the pressure does not stabilize within a few minutes, there could be air left in the pump. |
|---|--|

Note:

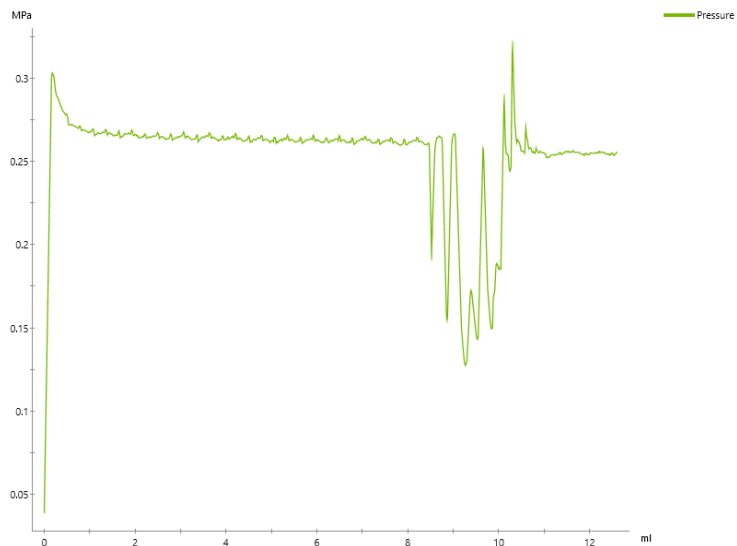
The pressure signal is considered stable if the fluctuation is no more than 5% up or down. See examples below.

Stable pressure (no air in the Pump):



Step	Action
------	--------

	Unstable pressure (air in the Pump)
--	-------------------------------------



- | | |
|---|--|
| 9 | If the pressure does not stabilize within a few minutes, repeat the procedures to prime the inlet tubing and purge the pump. If the pressure still does not stabilize, refer to the <i>ÄKTA go User Manual</i> for further instructions. |
|---|--|

Introduction

Details of the individual tests including purpose and required material are given in the method notes for each test method.

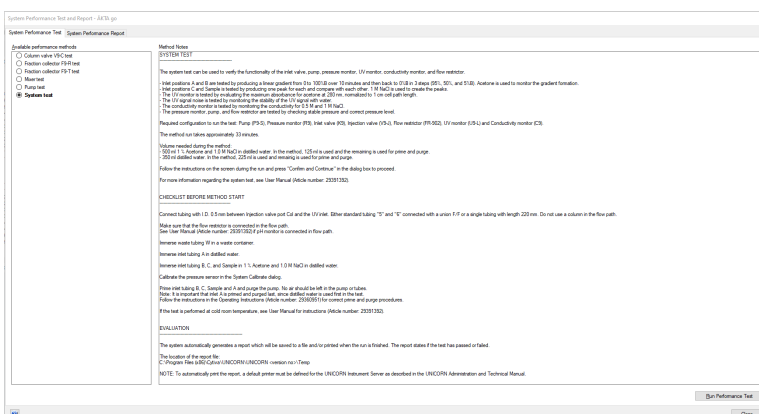
Note: The performance tests are always run without a column connected to the flow path.

Procedure

Performance tests are provided with the instrument configuration in the UNICORN software.

Follow the general steps below to run a performance test. Detailed requirements and procedures are shown on the screen when starting the test.

Step	Action
1	In the System Control module, select System → Performance Test and Report .
2	In the System Performance Test tab, in the list Available performance methods , select the test you want to run. Method notes for the selected test are shown in the right side panel.



Step	Action
------	--------

Note:

This dialog lists all tests that are available for the modules that can be installed in the ÅKTA go system. Attempting to run a test for a module that is not installed will generate an error message.

Note:

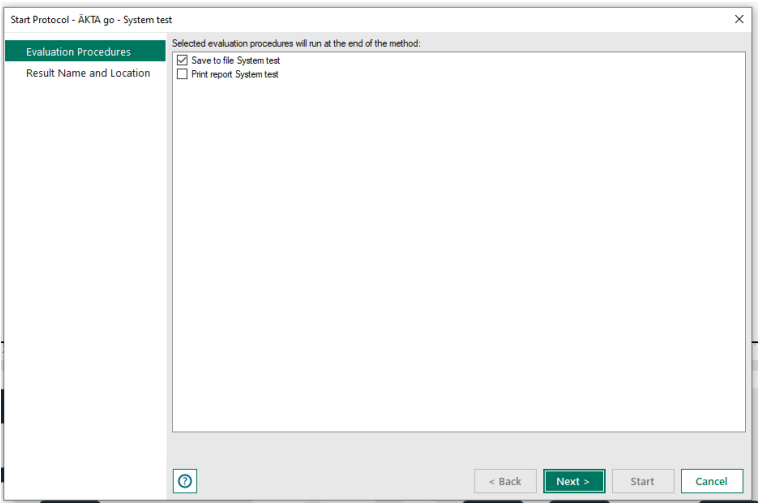
*The dialog for selecting and running a performance test includes a tab with the heading **System Performance Report**. This tab does not contain the reports for the performance tests.*

- 3 Read the information in the **Method Notes** panel carefully. Click **Run Performance Test**.

- 4 Tick the checkboxes **Save to file System test** and **Print report System test**. Click **Next >** to continue.

Note:

A printer must be installed on the UNICORN Instrument Server to be able to print the report (see the UNICORN Administration and Technical Manual for details).



- 5 Specify the details of the result file from the performance test. Click **Start** to start the performance test.

Note:

*The result file is separate from the test report. The report will be generated even if **No result** is selected.*

Step	Action
------	--------

- 6 Follow any instructions that are shown on the screen.
- 7 Check whether the test was passed or failed in the **System Performance Report**. The location of the report can be found in the **Method Notes** panel when selecting the test in the **System Performance Test** tab in the **System Control** module. The test result is stated at the top of the report. If the test was failed refer to *ÅKTA go User Manual* for possible causes.

Note: The progress of the performance test is shown in the **Chromatogram** pane in the **System Control** module.

5.4 Connect a column

Introduction

This section describes how to connect a column to the instrument, without introducing air into the flow path. Use a column holder to secure the column. Several types of column holders are available for ÄKTA go.



WARNING

To avoid exposing the column to excessive pressure, make sure that the pressure limit is set to the specified maximum pressure of the column. Before connecting a column to the ÄKTA go instrument, read the instructions for use of the column.



CAUTION

Cut injuries. The tubing cutter is very sharp and must be handled with care to avoid injuries.

Methods automatically include a pressure alarm based on the specifications of the chosen column. When performing a manual run, you must set the pressure limits yourself. See [Section 5.5 Pressure alarms, on page 78](#) for more information on pressure alarms.

Attach a column holder and connect a column

Follow the instructions below to connect a column to the instrument.

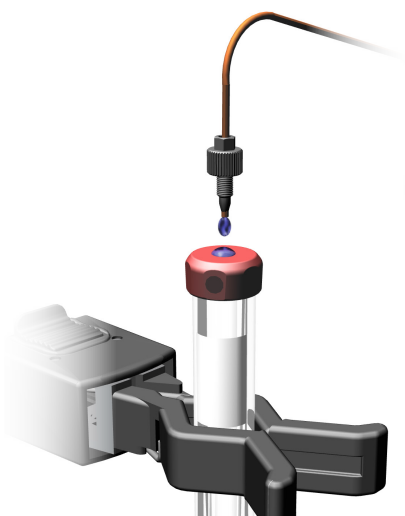
Note: *If the column is filled with 20% ethanol, do not use a salt-containing liquid when connecting it to the flow path because salt might precipitate in 20% ethanol.*

Step	Action
1	Attach an appropriate column holder to the rail on the instrument.
2	Attach the column to the column holder.
3	Cut a tubing of appropriate length to connect it between the injection valve and the top of the column. Connect this tubing to the column port Col of the injection valve.

Note:

Use red fingertights for columns with a red top and black for other columns.

- | Step | Action |
|------|--|
| 4 | <p>In the Process Picture pane, select Pump, enter a low System flow (e.g. 0.2 mL/min), and click Set.</p> <p><i>Result:</i></p> <p>A system flow of 0.2 mL/min starts.</p> <p>Note:</p> <p><i>If your system is equipped with a column valve, make sure to start a flow in the correct position in the column valve, and connect the column to those positions.</i></p> |
| 5 | <p>In the Process Picture pane, select Select limits in the Pressure pane. Choose an appropriate column or manually enter a pressure limit suitable for your column. Click Set.</p> <p><i>Result:</i></p> <p>A pressure limit suitable for your column is set.</p> |
| 6 | <p>Remove the stop plug from the top of the column.</p> |
| 7 | <p>When buffer is dripping from the tubing prepared in step 3, fill the top part of the column with buffer by letting it drip from the tube. When it is filled, remove the stop plug in the other end of the column and connect the tubing drop-to-drop to the top of the column.</p> |



Note:

Make sure that the connectors are properly tightened, but do not overtighten when connecting columns. Overtightening might rupture the connectors or squeeze the tubing and thereby result in high back pressure.

Step	Action
8	Cut a piece of tubing with appropriate length to connect the bottom of the column to the UV monitor. Remove the stop plug from the bottom of the column and connect this tubing in its place.
9	When buffer is dripping from the tubing connected to the bottom of the column, connect this piece of tubing to the UV monitor.

5.5 Pressure alarms

Introduction

The ÄKTA go instrument is equipped with a pressure monitor directly after the pump, which measures the highest pressure in the instrument. The advanced column valve, **V9-C**, adds two extra pressure sensors, one directly before the column and one directly after the column. These pressure monitors enable the measurement of pre-column pressure, which is the pressure on the column hardware, and delta-column pressure, which is the pressure on the packed bed.

Methods automatically include a pressure alarm based on the specifications of the chosen column. When performing a manual run, you must set the pressure limits yourself.

Note: *There is a default pressure limit of 2 MPa, but this limit does not protect all columns.*

Set a pressure alarm in a method

In a method, the pressure alarm is set in **Method Settings**. See [Section 6.1 Create a method, on page 81](#).

Set a pressure alarm in a manual run

In a manual run, the pressure alarm is set using the **Process Picture** pane or using the **Manual Instructions** box. In the **Process Picture**, click **Set limits** in the **Pressure** pane and enter a suitable pressure limit, or choose a column to get a suitable pressure limit. Click **Set**.

5.6 Prepare for a run at low temperature

Introduction

The viscosity of the liquids increase as the temperature decreases. Therefore, when using the instrument in a cold room or cold cabinet, decrease the flow rate and follow the precautions listed below.

Precautions concerning runs at cold room temperature



NOTICE

Avoid condensation. If ÄKTA go is kept in a cold room, cold cabinet or similar, keep it switched on in order to avoid condensation.



NOTICE

Avoid overheating. If ÄKTA go is kept in a cold cabinet and the cold cabinet is switched off, make sure to switch off ÄKTA go and keep the cold cabinet open to avoid overheating.



NOTICE

Place the computer in room temperature. If the product is placed in a cold room, use a cold room compatible computer or place the computer outside the cold room and use the Ethernet cable delivered with the instrument to connect to the computer.

Note: *When the instrument is kept in a cold room, it is important to retighten all tubing connectors and inlet connectors, because the plastics will shrink at a low temperature. Otherwise air might get into the flow path.*

Note: *Make sure that the instrument, buffers and sample have had time to reach the ambient temperature. When the instrument has reached the ambient temperature, calibrate all pressure sensors.*

6 Run a method

About this chapter

This chapter gives instructions on how to run your method.

In this chapter

Section	See page
6.1 Create a method	81
6.2 Prepare sample for loading	84
6.3 Start a method run	86
6.4 Monitor or interact with the run	87
6.5 Evaluate the run	88
6.6 Procedures after the run	90

6.1 Create a method

Introduction

This section describes how to create a new method in the **Method Editor**. There are more ways to create methods in UNICORN. Refer to the *UNICORN Method Manual* for more information.

Create a new method using the Method Editor

Follow the steps below to create a method using a predefined method template. The example given uses an **Affinity Chromatography** method. The **Method Settings** phase sets up parameters that are used throughout the method, such as unit for method base and flow rate.

- | Step | Action |
|------|--|
| 1 | Open the Method Editor module, click the New Method button and choose the system and a suitable predefined method. |
| 2 | Open the Phase Properties tab and confirm that all the selections made in the phases correspond to your intended method by following the steps below. |
| 3 | In the Method Settings phase, select column type to get suitable values, or enter correct values at Column Volume (column volume), Pressure Limit (pressure limit) and Flow Rate (flow rate). If using fraction collector F9-T select plate or tube type. |

The screenshot displays the 'Phase Properties' tab in the Method Editor. The left sidebar shows a vertical stack of phase buttons: 'Method Settings' (highlighted in blue), 'Equilibration', 'Sample Application', 'Column Wash - Wash Out Unbound Sample', 'Elution', and 'Re-Equilibration'. The main panel is divided into several sections:

- Method Settings:** Includes 'Show by Technique' (Affinity Chromatography (AC)), 'Column Type' (HiTrap HP, 1 ml), and a 'Column Type Properties...' button. Below this are 'Column Volume' (0.962 ml), 'Pressure Limit' (0.50 MPa), and a 'Use Flow Restrictor' checkbox.
- Flow Rate:** Includes a 'Flow Rate' input (1.00 ml/min) and checkboxes for 'Control to avoid Overpressure' and 'Reduce for Cold Room'.
- Fraction Collector:** Includes a 'Fraction Collector' dropdown (set to F9-T) and two 'Plate' dropdowns (both set to '96 deep well plate').
- Unit Selection:** Includes 'Method Base Unit' (CV) and 'Flow Rate Unit' (ml/min).
- Monitor Settings:** Includes a checkbox for 'Enable pH Monitoring'.
- Column Logbook:** Includes checkboxes for 'Enable logging of' and 'Cleaning in Place', and a 'Default Curve' dropdown (set to UV).

At the bottom of the interface are buttons for 'Delete', 'Save Phase...', and 'Duration & Variables'.

- | | |
|---|---|
| 4 | In the Equilibration phase, set a suitable volume and concentration of buffer B. |
|---|---|

Step	Action
------	--------

- | | |
|---|--|
| 5 | In the Sample Application phase, choose an appropriate sample application technique and volume. See Section 6.2 Prepare sample for loading, on page 84 for suitable sample application techniques for your sample volume. |
|---|--|

The screenshot displays the 'Sample Application' phase configuration in the ÅKTA go software. On the left, a vertical list of phases includes 'Method Settings', 'Equilibration', 'Sample Application' (which is highlighted in blue), 'Column Wash - Wash Out Unbound Sample', 'Elution', and 'Re-Equilibration'. Below this list are buttons for 'Delete', 'Save Phase...', and 'Duration & Variables'. The right pane is titled 'Sample Application' and contains several settings: 'Flow Rate' with a checkbox 'Keep the Flow Rate from the previous Phase' and a value of '1.00 ml/min'; 'Injection' with radio buttons for 'Loop Injection' and 'Pump Injection' (selected), and a 'Sample Inlet' dropdown set to 'Sample'; 'Inject fixed Sample Volume' set to '0.00 ml'; 'Finalize Sample Injection' set to '10.00 ml'; and 'Fractionation Settings' with radio buttons for 'Fraction Collector', 'Outlet Valve', 'Continue Ongoing', and 'No Fractionation' (selected). A note states: 'Select this option if no fractions are to be collected.'

Tip:

If you are using a sample loop, empty the loop with 3 to 5 volumes of buffer to make sure that all your sample fluid is flushed into the regular flow path.

Note:

For sample application using a Superloop™, refer to the ÅKTA go User Manual.

- | | |
|---|--|
| 6 | In the Column Wash phase, choose a suitable volume and concentration of buffer B, and whether a fraction collection is to be performed. |
| 7 | In the Elution phase, set how to elute the sample, the duration of the elution, the concentration of buffer B, and how the eluted sample will be collected. |
| 8 | In the Re-equilibration phase, choose a suitable volume and concentration of buffer B. |
| 9 | Click the save button, choose the system, name, and location for your method and click Save . |

Tip: *It is advisable to run a blank run, without sample, before running the method with sample. This makes sure that the column is clean and that the method and the system are set up properly.*

6.2 Prepare sample for loading

Introduction

This section describes how to prepare the sample for loading onto the column. This can be done using either a sample loop, a Superloop, or the pump. When using the pump to apply the sample, the sample inlet must first be primed.

Note: For sample application using a Superloop, refer to the ÄKTA go User Manual.

Sample application technique	Suitable volume
Sample loop	25 µL to 10 mL
Superloop	1 to 150 mL
Pump	From 5 mL

Prepare sample for application using a sample loop

Follow the steps below to prepare for sample application using a sample loop.

Step	Action
1	Connect a suitable sample loop to the injection valve ports LoopF (fill) and LoopE (empty).
2	Fill a syringe with sample.
3	Connect the syringe to the injection valve port Syr .
4	In the Process Picture , make sure that the injection valve is in position Load .

Note:

Load is the default position for the valve.

5	Load sample into the sample loop. To avoid sample loss due to siphoning, leave the syringe in the port until the sample has been injected onto the column during the run.
---	---

Tip:

When repeatability is important, overload the loop with 3 to 5 volumes of sample to make sure that the loop is completely filled.

Note:

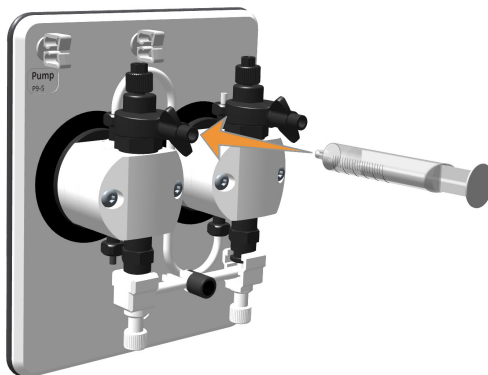
Make sure the waste bottle is located below the injection valve, to avoid waste going back into the sample loop.

Prepare sample for application using the pump

Follow the steps below to prepare for sample application using the pump.

Step	Action
------	--------

- | | |
|---|--|
| 1 | Make sure that the sample inlet tubing that is to be used during the run is placed in the sample container. |
| 2 | In the Process Picture pane, click on Inlet Valve and select Sample .
<i>Result:</i>
The inlet valve opens the Sample inlet. |
| 3 | Connect a 25 to 30 mL syringe to the purge valve of one of the pump heads. |



- | | |
|---|--|
| 4 | Open the purge valve by turning it counter-clockwise one and a half turns. Draw liquid slowly into the 25 to 30 mL syringe until the liquid reaches the inlet valve. |
| 5 | Close the purge valve by turning it clockwise. Disconnect the syringe and discard its contents. |

6.3 Start a method run

Introduction

This section describes how to start a run using a previously created method. For further information on method creation, please refer to *UNICORN Method Manual*.



Prerequisites

Make sure that the system is correctly prepared. Confirm the following:

- The sample inlet is primed or the sample loop is loaded.
- The column is properly connected.
- The pressure alarm has an appropriate limit set.
- That there is no air in the system.
- The buffer inlet tubing(s) is immersed in correct buffer vessels.
- All waste tubing is immersed in appropriate waste vessels that have sufficient empty volume.
- No tubing is twisted and the flow path is free from leakage.
- If a fraction collector is being used, make sure that the correct plates and/or tubes are used and loaded.

Start a method run

Follow the steps below to start the method run.

Step	Action
1	Open the System Control module and click the button Open Method Navigator . 
2	Select the method to run, and click the Run button. 
3	Step through the displayed pages in the Start Protocol , add requested input and make appropriate changes if necessary. Click Next .
4	Click Start on the last page of the Start Protocol .

6.5 Evaluate the run

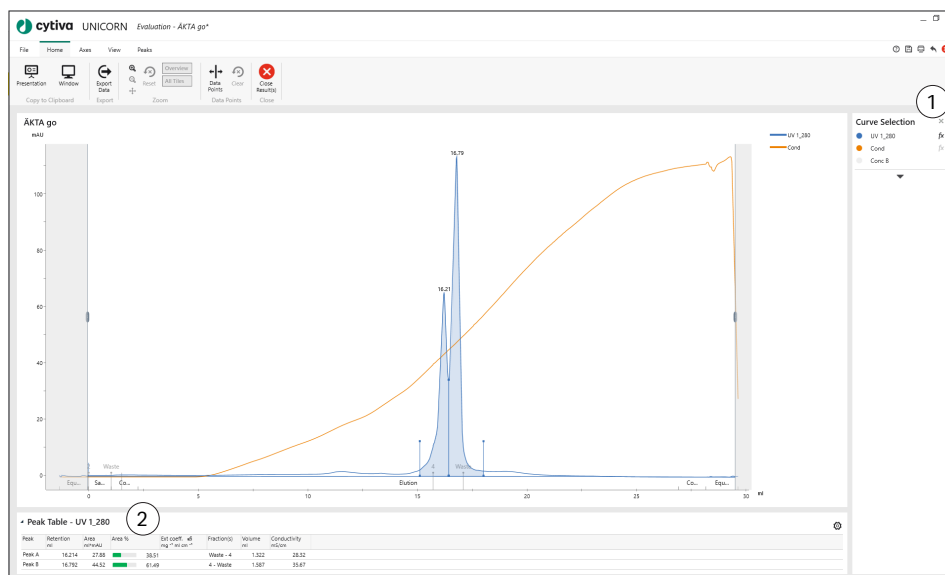
Introduction

After the run is completed, the results can be evaluated using the **Evaluation** module. To open the results, use the following steps. For more information on evaluating results, refer to the UNICORN user documentation.

Step	Action
1	In the Evaluation module, click Result . Browse or search for your result file.
2	Check in the preview that you have the correct result file and click Open .
3	To save an evaluation file, click File → Save .

Illustration of an evaluated run

The image below shows an example of the features available in the **Evaluation** module in UNICORN.



Part	Function
1	To integrate or remove integration click the fx button next to the curve name. To adjust integration parameters go to the Peak tab.

Part	Function
2	Data from the integration is shown in a Peak table below the chromatogram. Customize the Peak data table by clicking the settings icon on the top right of the Peak table.

6.6 Procedures after the run

Introduction

After the run, the instrument and column should be cleaned to prevent bacterial growth, sample contamination in the next run, and column clogging.

This section describes how to clean the column using a **Column CIP** (Cleaning-In-Place) method, prepare the column and the instrument for storage, and how to shut down the system.

Create a CIP method

To create a **Column CIP** method, use the following procedure.

Step	Action
1	Open the Method Editor module and click the Create new method button.
2	Select Column CIP method .
3	In the method settings, select the column being used.
4	In Column CIP phase, click get suggested steps .
	Note: <i>The suggested steps are only available for some techniques and might not be optimized for your column. Read the instructions provided with the column to get the recommended CIP steps.</i>
5	Click the save button, choose the system, name, and location for your method and click Save .

Run a column CIP

To clean the column using the **Column CIP** method, use the following procedure.

Step	Action
1	Open the System Control module and click the Open Method Navigator button.
2	Select the Column CIP method created above and click the Run button.
3	Make sure that the inlet tubings are in the correct solutions.
4	Click Start .

Column storage

If the column is not going to be used for a couple of days or longer, it must be placed in a storage solution (e.g. 20% ethanol) after a **Column CIP** run. Refer to the instructions for your column for specific storage instructions. To place the column in a storage solution, use the following procedure.

Note: *If the column is in a salt-containing buffer, equilibrate the column in water before putting it in 20% ethanol, otherwise salt might precipitate in the column.*

Step	Action
1	Place the inlet tubing in the storage solution (20% ethanol).
2	In the Process Picture , click on Pump .
3	Enter a flow rate of half the recommended flow rate of your column.
4	Click on Timer and select Volume .
5	Under Volume , enter 4 times the column volume. Click Start .
6	When the run has finished, remove the column, plug the top and bottom of the column, and place in a refrigerator.

System storage

If the system is not going to be used for a couple of days or longer, clean the system using the **System CIP** method (see [Section 7.2 Perform system Cleaning-In-Place \(CIP\), on page 95](#)) and then put the system in storage solution. Start by removing the cleaned column from the system, see above. To put the system in storage, follow the steps below.

Step	Action
1	Clean the system using a System CIP method created for the system, see Section 7.2 Perform system Cleaning-In-Place (CIP), on page 95 .
2	Place all inlet tubing in storage solution (20% ethanol or 2% benzyl alcohol).
3	In the Process Picture , click on Pump .
4	Tick the Wash box for all inlet tubing in storage solution and click Pump Wash .
5	Enter a flow rate of 10 mL/min.
6	Click on Timer , select Volume , and enter 25 mL.
7	Decrease the flow rate to 2 mL/min and run a fractionation for two minutes to fill the fraction collector tubing with storage solution.

Step	Action
8	Replace the rinsing solution with 20% ethanol. See Section 4.2.3 Prepare the pump rinsing system, on page 50 .
Note:	<i>If the system is equipped with a column valve, make sure all positions are cleaned with storage solution.</i>
Note:	<i>Make sure to clean all applicable positions and ports of the outlet valve.</i>

Shut down the system

Follow the steps below to shut down the system.

Step	Action
1	Select Exit UNICORN from the File menu in any module in the UNICORN software.
2	Press the On/Off button on the instrument control panel.



WARNING

Power is still supplied to some internal electronics circuits when the instrument is switched off using the On/Off button. Disconnect the instrument from the power supply before maintenance or service.



NOTICE

Damage to RTU. To shut down the system, make sure to press the On/Off button for a short time. When the button is pressed for an extended period of time, the system promptly shut down, which may damage the Real-Time Unit (RTU).

7 Maintenance

About this chapter

This chapter provides information on how to perform basic maintenance procedures. For a complete list of maintenance procedures, refer to the *ÄKTA go User Manual*.



WARNING

All maintenance procedures inside the instrument chassis must be performed by a Cytiva service representative.

In this chapter

Section		See page
7.1	Clean the instrument externally	94
7.2	Perform system Cleaning-In-Place (CIP)	95
7.3	Replace pump rinsing liquid	99
7.4	Replace the main fuses	100

7.1 Clean the instrument externally



CAUTION

Disconnect power. Always disconnect power from the instrument before performing any maintenance task.

Maintenance interval

Clean the instrument externally when required.

Required material

The following materials are required:

- Cloth
- Mild cleaning agent or 20% ethanol

Instruction

Follow the steps below to clean the instrument externally.

Step	Action
1	Turn off the instrument and disconnect the power cord from the wall socket.
2	Wipe the surface with a damp cloth. Wipe off stains using a mild cleaning agent or 20% ethanol. Wipe off any excess.
3	Remove the top and bottom tray for cleaning. Put them back once they are cleaned.
4	Let the instrument dry completely before using it.

7.2 Perform system Cleaning-In-Place (CIP)

Introduction

The **System CIP** method is used to clean the flow path.



WARNING

Explosion hazard due to leakage of flammable liquid. Make sure there is no leakage in the flow path of the instrument during the cleaning with a flammable liquid, such as ethanol. Carefully inspect the flow path for leakage, including the waste tubing and tighten any connectors if necessary.



WARNING

Hazardous biological agents during run. When using hazardous biological agents, perform a system cleaning to flush the entire system tubing with storage solution (e.g. NaOH) followed by a neutral buffer and finally distilled water, before service and maintenance.



CAUTION

Hazardous substances. When using hazardous chemical and biological agents, take all suitable protective measures, such as wearing protective glasses and gloves resistant to the substances used. Follow local and/or national regulations for safe operation, maintenance and decommissioning of the equipment.

Tip: *If hazardous chemicals are used for system or column cleaning, wash the system or columns with a neutral solution in the last phase or step of the system CIP method.*

Maintenance interval

Perform a system CIP when required, for example between runs where different samples are used. This is important to prevent cross-contamination and bacterial growth in the instrument.

Required material

The following materials are required:

- Appropriate cleaning solutions (1 M NaOH, 0.1 M NaCl, buffer solution, and distilled water).

- Syringe, 25 to 30 mL

Create a System CIP method

The default **System CIP** method comprises of phases that cleans the system with water, 1 M NaOH, 0.1 M NaCl, and buffer. Default inlets are **B**, **C**, **Sample**, and **A**, respectively. The inlets used can be changed in each phase. Follow the steps below to create a **System CIP** method.

Note: *The pH valve and the pH electrode are not included in the **System CIP** method. For information about cleaning of the pH valve and the pH electrode, refer to the ÄKTA go User Manual.*

Step	Action
------	--------

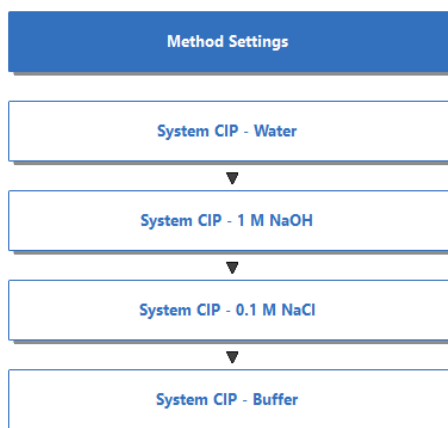
- | | |
|---|--|
| 1 | In the Method Editor module, click the Create New Method button. |
|---|--|



- | | |
|---|---|
| 2 | In the New Method dialog, select System and in the Predefined Method drop-down list select System CIP . Click OK . |
|---|---|

Result:

A **System CIP** method is opened, comprising the **Method Settings** phase and four **System CIP** phases. Each **System CIP** phase uses one cleaning solution.



- | | |
|---|---|
| 3 | If applicable, add additional System CIP phases to the method by dragging and dropping them from the Phase Library . |
| 4 | Click the Save button. |
| 5 | In the Save As dialog, select a target folder, enter the Name for the method, select a System from the list, and click Save . |

Perform a System CIP

Follow the steps below to run a **System CIP** method.

Step	Action
1	Remove the column tubing and connect tubing between the Col port on the injection valve and the UV monitor. Note: <i>If you have a column valve, connect bypass tubing to all column positions.</i>
2	Prepare cleaning solutions and immerse the selected inlet tubing in the solutions. Note: <i>The default solutions to use are buffer for inlet A, water for inlet B, 1 M NaOH for inlet C, and 0.1 M NaCl for the Sample inlet.</i>
3	In the System Control module, select the System CIP method created above, and start the run.
4	For complete cleaning of the flow path, clean the manual injection port of the injection valve manually, see the instructions below.

Clean the syringe port of the injection valve

Follow the steps below to manually clean the syringe port and the sample loop.

Step	Action
1	In the Process Picture make sure that the injection valve is in position Load . Note: <i>Load is the default position for the valve.</i>
2	Connect a suitable sample loop to the injection valve ports LoopF (fill) and LoopE (empty). Note: <i>Do not use a Superloop when cleaning the Injection valve.</i>
3	Make sure there is waste tubing connected to the injection valve port W2 and W1 and that they are secured in a waste container.

Step	Action
------	--------

- | | |
|---|---|
| 4 | Fill a syringe with approximately 10 mL of an appropriate cleaning solution (1 M NaOH or buffer solution). Connect the syringe to injection valve port Syr , and inject the cleaning solution. |
|---|---|



- | | |
|---|---|
| 5 | Fill a syringe with distilled water. Connect the syringe to injection valve port Syr , and inject the distilled water. |
|---|---|

7.3 Replace pump rinsing liquid

Replace pump rinsing liquid once a week if you are using 20% ethanol and daily if you are using an aqueous buffer. For instructions, see [Section 4.2.3 Prepare the pump rinsing system, on page 50](#).

7.4 Replace the main fuses



WARNING

Disconnect power. Always disconnect power from the instrument before replacing fuses.



WARNING

For continued protection from fire hazard, replace only with same type and rating of fuse.



WARNING

If a fuse requires repeated replacement, do not continue to use the instrument. Contact an authorized service engineer.

Required tools

Tool	Dimension
Flat screwdriver	2 to 3 mm

Procedure

Follow the instructions below to replace the fuses.

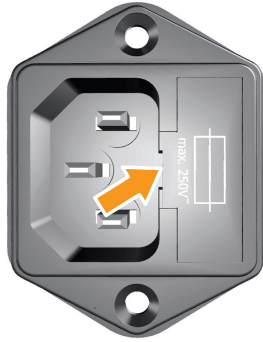
Note: *There are two identical fuses at the mains power inlet. Replace both fuses even if only one is blown.*

Step	Action
------	--------

- | | |
|---|---|
| 1 | Turn off the instrument. |
| 2 | Disconnect the power cord from the power inlet. |

Step	Action
------	--------

- | | |
|---|---|
| 3 | Use a small flat-bladed screwdriver to prise open the fuse holder cover on the power inlet. |
|---|---|



- | | |
|---|--|
| 4 | Pull the fuse holder out of the mains connector panel by hand. |
| 5 | Remove the fuses from the fuse holder. |
| 6 | Fit new fuses with size 5 x 20 mm and rating T4AL 250 V. |
| 7 | Replace the fuse holder in the power inlet. |



NOTICE

When replacing fuses, make sure that the fuse holder is pushed fully into position.

- | | |
|---|---------------------------|
| 8 | Reconnect the power cord. |
|---|---------------------------|

8 Troubleshooting

About this chapter

This chapter provides a list of the most commonly encountered problems that might occur when operating ÄKTA go. For a more comprehensive list and a more detailed description of the actions to take, refer to *ÄKTA go User Manual*.

Problem	Possible cause	Corrective action
Spike in the UV signal	The Flow restrictor has been removed after the UV cell	Replace the Flow restrictor back in the flow path after the Conductivity monitor
	There is air in the system	Prime the inlets and purge the pump, see Section 5.2 Prime inlets and purge pump heads, on page 66
Unstable conductivity signal	There is air in the pump	Prime the inlets and purge the pump, see Section 5.2 Prime inlets and purge pump heads, on page 66
Unstable pressure	There is air in the pump	Prime the inlets and purge the pump, see Section 5.2 Prime inlets and purge pump heads, on page 66
Pressure alarm issued	The pressure is too high, possibly due to running the system at a flow rate that is too high for the column used	Lower the flow rate
	There is a kink in the tubing, overtightened tubing connections, or precipitation in the flow path	Replace the tubing. If replacement of tubing does not fix the problem, refer to the <i>ÄKTA go User Manual</i> , to continue troubleshooting

Problem	Possible cause	Corrective action
Difficulty connecting to the system	IP addresses are on different subnets	Make sure the instrument IP address and the computer IP address are on the same subnet, refer to <i>UNICORN Administration and Technical manual</i> . The instrument IP address can be seen on the instrument control panel when the instrument is powered on but does not have connection to UNICORN.
	UNICORN instrument server is not started	Restart the computer
The text in UNICORN is large and cannot be read because it is truncated	Scaling in the Windows operating system has been changed	Set the zoom level in the Windows operating system to 100% and restart the computer

System error report

When you request troubleshooting assistance from Cytiva, you should generate a **System error report** (system error report) and submit it to your service representative.

Follow the instructions below to generate a **System error report**.

Step	Action
1	Select System → Create System Error Report in the System Control module. Users with sufficient access rights can also create a System error report from the Administration module.
2	Step through the report wizard using the Next and Back buttons. Provide information as requested at each step. Add results, methods and logs as appropriate.
3	Save the report in the default folder. The report is saved as a zip file with the name Report_YYYYMMDDnn.zip .
4	Submit the file to your Cytiva service representative.
	Note: <i>The file may be large (> 15 Mb).</i>

9 Reference information

About this chapter

This chapter contains reference information.

In this chapter

Section		See page
9.1	System specifications	105
9.2	Chemical resistance specifications	106
9.3	Recycling information	108
9.4	Regulatory information	109

9.1 System specifications

Technical specifications

Parameter	Specification
System configuration	Benchtop system, external computer
Flow rate range	0.01 to 25 mL/min
Pressure range	0 to 5 MPa (0 to 50 bar)
Control system	ÄKTA go 1.0 UNICORN 7.4 or later version ÄKTA go 2.0 UNICORN 7.6 or later version ÄKTA go 3.0 UNICORN 7.6 or later version
Connection between PC and instrument	Ethernet
Dimensions (width × height × depth)	335 × 482 × 464 mm (depth without tray 451 mm, depth without modules 380 mm)
Weight (excluding computer, columns, buffer bottles)	27 kg
Supply voltage	100 to 240 V~ autorange
Maximum voltage fluctuation	± 10% from the nominal voltage
Frequency	50/60 Hz
Power consumption	Rated max 300 VA ¹ Max with all options 150 W ² Typical 100 W Power-save < 20 W
Fuse rating	T4AL 250 V
Enclosure protective class	IP 21
Acoustic noise level	< 60 dB(A)

¹ ÄKTA go can deliver 300 VA.

² ÄKTA go equipped with all options consume 150 W.

Battery information

The instrument version with Real-Time Unit contains a lithium backup battery. The battery cannot be replaced by the user.

9.2 Chemical resistance specifications

Introduction

This section provides detailed information about chemical resistance of the ÄKTA go instrument to some common aggressive chemicals used in liquid chromatography. For information regarding chemicals not listed in this section, contact your Cytiva representative.

Note: Refer to Safety Data Sheets (SDS) for information regarding characteristics, human and environmental risks and preventive measures for chemicals used. Make sure that you have the SDS available from your chemical distributor and/or databases on the internet.

Considerations

The information in this section applies to the ÄKTA go flow path. After using organic solvents, restore the system in distilled water (e.g. overnight) or storage solution (20% ethanol or 2% benzyl alcohol) as soon as possible. To spare flow path material, do not leave it exposed to these chemicals more than required for the chromatographic runs. The definition for long-term use is approximately 8 hours/day, 5 days/week. Short-term use is defined as less than 2 hours.

Recommended solutions for the pump rinsing system are aqueous buffers or 20% ethanol. Do not use other chemicals in the pump rinsing system.

Chemical resistance, long-term use

The following chemicals are suitable for continuous (long-term) use.

Chemical	Concentration	CAS no	EC no
Aqueous buffers, pH 2 to 12	N/A	N/A	N/A
Acetone	10%	67-64-1	200-662-2
Acetonitrile	83%	75-05-8	200-835-2
Ammonia	30%	7664-41-7	231-635-3
Ammonium chloride	2 M	12125-02-9	235-186-4
Ammonium sulfate	3 M	7783-20-2	231-984-1
Arginine	2 M	74-79-3	200-811-1
Benzyl alcohol ¹	2%	100-51-6	202-859-9
Dimethyl sulfoxide (DMSO)	5%	67-68-5	200-664-3
Dithioerythritol (DTE)	100 mM	6892-68-8	229-998-8
Dithiothreitol (DTT)	100 mM	3483-12-3	222-468-7

Chemical	Concentration	CAS no	EC no
Ethanol ¹	20%	75-08-1	200-837-3
Ethylene glycol	50%	107-21-1	203-473-3
Ethylenediaminetetraacetic acid (EDTA)	100 mM	60-00-4	200-449-4
Glycerol	50%	56-81-5	200-289-5
Guanidinium hydrochloride	6 M	50-01-1	200-002-3
Mercaptoethanol	20 mM	37482-11-4	253-523-3
Phosphoric acid	0.1 M	7664-38-2	231-633-2
Potassium chloride	4 M	7447-40-7	231-211-8
Sodium dodecyl sulfate (SDS)	1%	151-21-3	205-788-1
Sodium hydroxide	0.01 M	1310-73-2	215-185-5
Tween 20	1%	9005-64-5	500-018-3
Urea	8 M	57-13-6	200-315-5

¹ Suitable storage solution

Chemical resistance, short-term use

The following chemicals are suitable for up to 2 h contact time at room temperature.

Chemical	Concentration	CAS no	EC no
Acetic acid	70%	75-05-8	200-835-2
Decon 90	10%	N/A	N/A
Ethanol	100%	75-08-1	200-837-3
Hydrochloric acid	0.1 M	7647-01-0	231-595-7
Isopropanol	100%	67-63-0	200-661-7
Methanol	100%	67-56-1	200-659-6
Sodium hydroxide	2 M	1310-73-2	215-185-5
Sodium hydroxide/ethanol	1 M/40%	N/A	N/A
Sodium chloride	4 M	7647-14-5	231-598-3
Sodium hypochlorite	10%	7681-52-9	231-668-3

9.3 Recycling information

Introduction

This section contains information about the decommissioning of the product.



CAUTION

Always use appropriate personal protective equipment when decommissioning the equipment.

Decontamination

The product must be decontaminated before decommissioning. All local regulations must be followed with regard to scrapping of the equipment.

Disposal of the product

When taking the product out of service, the different materials must be separated and recycled according to national and local environmental regulations.

Recycling of hazardous substances

The product contains hazardous substances. Detailed information is available from your Cytiva representative.

Disposal of electrical components



Waste electrical and electronic equipment must not be disposed of as unsorted municipal waste and must be collected separately. Contact an authorized representative of the manufacturer for information concerning the decommissioning of the equipment.

Disposal of batteries

Waste batteries and accumulators must not be disposed of as unsorted municipal waste and must be collected separately. Follow applicable local regulations for recycling of batteries and accumulators.

9.4 Regulatory information

Introduction

This section lists the regulations and standards that apply to the product. Your system is marked or listed according to the applicable regulatory requirements for your region. Local language translations are only provided according to regulatory requirements.

In this section

Section		See page
9.4.1	Contact information	110
9.4.2	European Union and European Economic Area	111
9.4.3	Great Britain	112
9.4.4	Eurasian Economic Union (Евразийский экономический союз)	113
9.4.5	North America	115
9.4.6	General regulatory statements	116
9.4.7	South Korea	117
9.4.8	China	118
9.4.9	Health and Safety Declaration Form	121

9.4.1 Contact information

Contact information for support

To find local contact information for support and sending troubleshooting reports, visit cytiva.com/contact.

Manufacturing information

The table below summarizes the required manufacturing information.

Requirement	Information
Name and address of manufacturer	Cytiva Sweden AB Björkgatan 30 SE 751 84 Uppsala Sweden
Telephone number of manufacturer	+ 46 771 400 600

9.4.2 European Union and European Economic Area

Introduction

This section describes regulatory information for the European Union and European Economic Area that applies to the product.

Conformity with EU Directives

See the EU Declaration of Conformity for the directives and regulations that apply for the CE marking.

If not included with the product, a copy of the EU Declaration of Conformity is available on request.

CE marking



The CE marking and the corresponding EU Declaration of Conformity is valid for the product when it is:

- used according to the *Operating Instructions* or user manuals, and
- used in the same state as it was delivered, except for alterations described in the *Operating Instructions* or user manuals.

9.4.3 Great Britain

Introduction

This section describes regulatory information for Great Britain that applies to the product.

Conformity with UK Regulations

See the UK Declaration of Conformity for the regulations that apply for the UKCA marking.

If not included with the product, a copy of the UK Declaration of Conformity is available on request.

UKCA marking



The UKCA marking and the corresponding UK Declaration of Conformity is valid for the product when it is:

- used according to the *Operating Instructions* or user manuals, and
- used in the same state as it was delivered, except for alterations described in the *Operating Instructions* or user manuals.

9.4.4 Eurasian Economic Union (Евразийский экономический союз)

This section describes the information that applies to the product in the Eurasian Economic Union (the Russian Federation, the Republic of Armenia, the Republic of Belarus, the Republic of Kazakhstan, and the Kyrgyz Republic).

Introduction

This section provides information in accordance with the requirements of the Technical Regulations of the Customs Union and (or) the Eurasian Economic Union.

Введение

В данном разделе приведена информация согласно требованиям Технических регламентов Таможенного союза и (или) Евразийского экономического союза.

Manufacturer and importer information

The following table provides summary information about the manufacturer and importer, in accordance with the requirements of the Technical Regulations of the Customs Union and (or) the Eurasian Economic Union.

Requirement	Information
Name, address and telephone number of manufacturer	See <i>Manufacturing information</i>
Importer and/or company for obtaining information about importer	<p>Cytiva RUS LLC</p> <p>109004, Moscow</p> <p>internal city area Tagansky municipal district</p> <p>Stanislavsky str., 21, building 5, premises I, offices 24,25,29</p> <p>Russian Federation</p> <p>Telephone: +7 985 192 75 37</p> <p>E-mail: rucis@cytiva.com</p>

Информация о производителе и импортере

В следующей таблице приводится сводная информация о производителе и импортере, согласно требованиям Технических регламентов Таможенного союза и (или) Евразийского экономического союза.

Требование	Информация
Наименование, адрес и номер телефона производителя	См. Информацию об изготовлении
Импортёр и/или лицо для получения информации об импортёре	<p>ООО "Цитива РУС"</p> <p>109004, г. Москва</p> <p>вн. тер. г. муниципальный округ Таганский</p> <p>ул. Станиславского, д. 21 стр. 5, помещ. I, ком. 24,25,29</p> <p>Российская Федерация</p> <p>Телефон: +7 985 192 75 37</p> <p>Адрес электронной почты: rucis@cytiva.com</p>

Description of symbol on the nameplate

Описание символов на заводской табличке



This Eurasian compliance mark indicates that the product is approved for use on the markets of the Member States of the Customs Union of the Eurasian Economic Union

Данный знак о Евразийском соответствии указывает, что изделие одобрено для использования на рынках государств-членов Таможенного союза Евразийского экономического союза

9.4.5 North America

Introduction

This section describes the information that applies to the product in the USA and Canada.

FCC compliance

This device complies with part 15 of the FCC Rules. Operation is subject to the following two conditions: (1) This device may not cause harmful interference, and (2) this device must accept any interference received, including interference that may cause undesired operation.

Note: *The user is cautioned that any changes or modifications not expressly approved by Cytiva could void the user's authority to operate the equipment.*

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

9.4.6 General regulatory statements

Introduction

This section shows regulatory statements that are applicable to more than one geographical region.

EMC emission, CISPR 11: Group 1, Class A statement



NOTICE

This equipment is not intended for use in residential environments and may not provide adequate protection to radio reception in such environments.

9.4.7 South Korea

Introduction

This section describes the regulatory information to comply with the Korean technical regulations.

Compliance statement

**NOTICE**

Class A equipment (equipment for business use).

This equipment has been evaluated for its suitability for use in a business environment.

When used in a residential environment, there is a concern of radio interference.

**유의사항**

A급 기기(업무용 방송통신기자재)

이 기기는 업무용 환경에서 사용할 목적으로 적합성평가를 받은 기기

로서 가정용 환경에서 사용하는 경우 전파간섭의 우려가 있습니다.

9.4.8 China

This section describes the information that applies to the product in China.

有害物质声明 (DoHS)

Declaration of Hazardous Substances

(DoHS)

根据 SJ/T11364-2014 《电子电气产品有害物质限制使用标识要求》特提供如下有关污染控制方面的信息。

The following product pollution control information is provided according to SJ/T11364-2014 Marking for Restriction of Hazardous Substances caused by electrical and electronic products.

电子信息产品污染控制标志说明

Explanation of Pollution Control Label



该标志表明本产品含有超过中国标准 GB/T 26572 《电子电气产品中限用物质的限量要求》中限量的有害物质。标志中的数字为本产品的环保使用期，表明本产品在不正常使用的条件下，有毒有害物质不会发生外泄或突变，用户使用本产品不会对环境造成严重污染或对其人身、财产造成严重损害的期限。单位为年。

为保证所声明的环保使用期限，应按产品手册中所规定的环境条件和方法进行正常使用，并严格遵守产品维修手册中规定的定期维修和保养要求。

产品中的消耗件和某些零部件可能有其单独的环保使用期限标志，并且其环保使用期限有可能比整个产品本身的环保使用期限短。应到期按产品维修程序更换那些消耗件和零部件，以保证所声明的整个产品的环保使用期限。

本产品在使用寿命结束时不可作为普通生活垃圾处理，应被单独收集妥善处理。

This symbol indicates the product contains hazardous materials in excess of the limits established by the Chinese standard GB/T 26572 Requirements of concentration limits for certain restricted substances in electrical and electronic products. The number in the symbol is the Environment-friendly Use Period (EFUP), which indicates the period during which the hazardous substances contained in electrical and electronic products will not leak or mutate under normal operating conditions so that the use of such electrical and electronic products will not result in any severe environmental pollution, any bodily injury or damage to any assets. The unit of the period is "Year".

In order to maintain the declared EFUP, the product shall be operated normally according to the instructions and environmental conditions as defined in the product manual, and periodic maintenance schedules specified in Product Maintenance Procedures shall be followed strictly.

Consumables or certain parts may have their own label with an EFUP value less than the product. Periodic replacement of those consumables or parts to maintain the declared EFUP shall be done in accordance with the Product Maintenance Procedures.

This product must not be disposed of as unsorted municipal waste, and must be collected separately and handled properly after decommissioning.

有害物质的名称及含量
Name and Concentration of
Hazardous Substances

产品中有害物质的名称及含量
Table of Hazardous Substances' Name and Concentration

部件名称 Component name	有害物质 Hazardous substance					
	铅 (Pb)	汞 (Hg)	镉 (Cd)	六价铬 (Cr(VI))	多溴联苯 (PBB)	多溴二苯醚 (PBDE)
29375260	X	0	0	0	0	0

- 0:** 表示该有害物质在该部件所有均质材料中的含量均在 GB/T 26572 规定的限量要求以下。

X: 表示该有害物质至少在该部件的某一均质材料中的含量超出 GB/T 26572 规定的限量要求。

 - 此表所列数据为发布时所能获得的最佳信息。
- 0:** Indicates that this hazardous substance contained in all of the homogeneous materials for this part is below the limit requirement in GB/T 26572.

X: Indicates that this hazardous substance contained in at least one of the homogeneous materials used for this part is above the limit requirement in GB/T 26572

 - Data listed in the table represents best information available at the time of publication.

9.4.9 Health and Safety Declaration Form

On site service



On Site Service Health & Safety Declaration Form

Service Ticket #:	
--------------------------	--

To make the mutual protection and safety of Cytiva service personnel and our customers, all equipment and work areas must be clean and free of any hazardous contaminants before a Service Engineer starts a repair. To avoid delays in the servicing of your equipment, complete this checklist and present it to the Service Engineer upon arrival. Equipment and/or work areas not sufficiently cleaned, accessible and safe for an engineer may lead to delays in servicing the equipment and could be subject to additional charges.

Yes	No	Review the actions below and answer "Yes" or "No". Provide explanation for any "No" answers in box below.
<input type="radio"/>	<input type="radio"/>	Instrument has been cleaned of hazardous substances. Rinse tubing or piping, wipe down scanner surfaces, or otherwise make sure removal of any dangerous residue. Make sure the area around the instrument is clean. If radioactivity has been used, perform a wipe test or other suitable survey.
<input type="radio"/>	<input type="radio"/>	Adequate space and clearance is provided to allow safe access for instrument service, repair or installation. In some cases this may require customer to move equipment from normal operating location prior to Cytiva arrival.
<input type="radio"/>	<input type="radio"/>	Consumables, such as columns or gels, have been removed or isolated from the instrument and from any area that may impede access to the instrument.
<input type="radio"/>	<input type="radio"/>	All buffer / waste vessels are labeled. Excess containers have been removed from the area to provide access.
Provide explanation for any "No" answers here:		
Equipment type / Product No:		Serial No:
I hereby confirm that the equipment specified above has been cleaned to remove any hazardous substances and that the area has been made safe and accessible.		
Name:		Company or institution:
Position or job title:		Date (YYYY/MM/DD):
Signed:		

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

© 2020 Cytiva.

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact).
 28980026 AD 04/2020

Product return or servicing



Health & Safety Declaration Form for Product Return or Servicing

Return authorization number:		and/or Service Ticket/Request:	
-------------------------------------	--	---------------------------------------	--

To make sure the mutual protection and safety of Cytiva personnel, our customers, transportation personnel and our environment, all equipment must be clean and free of any hazardous contaminants before shipping to Cytiva. To avoid delays in the processing of your equipment, complete this checklist and include it with your return.

1. Note that items will NOT be accepted for servicing or return without this form
2. Equipment which is not sufficiently cleaned prior to return to Cytiva may lead to delays in servicing the equipment and could be subject to additional charges
3. Visible contamination will be assumed hazardous and additional cleaning and decontamination charges will be applied

Yes	No	Specify if the equipment has been in contact with any of the following:	
<input type="radio"/>	<input type="radio"/>	Radioactivity (specify)	
<input type="radio"/>	<input type="radio"/>	Infectious or hazardous biological substances (specify)	
<input type="radio"/>	<input type="radio"/>	Other Hazardous Chemicals (specify)	
Equipment must be decontaminated prior to service / return. Provide a telephone number where Cytiva can contact you for additional information concerning the system / equipment.			
Telephone No:			
Liquid and/or gas in equipment is:		<input type="checkbox"/>	Water
		<input type="checkbox"/>	Ethanol
		<input type="checkbox"/>	None, empty
		<input type="checkbox"/>	Argon, Helium, Nitrogen
		<input type="checkbox"/>	Liquid Nitrogen
		<input type="checkbox"/>	Other, specify
Equipment type / Product No:		Serial No:	
I hereby confirm that the equipment specified above has been cleaned to remove any hazardous substances and that the area has been made safe and accessible.			
Name:		Company or Institution:	
Position or job title:		Date (YYYY/MM/DD)	
Signed:			

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

© 2020 Cytiva.

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact).
29980027 AD 04/2020

To receive a return authorization number or service number, call local technical support or customer service.

Index

A

ÄKTA go system, 105
 specifications, 105
 Ambient environment, 39

C

CE, 111
 conformity, 111
 marking, 111
 CIP, 95
 system CIP, 95
 Connect system units, 45
 Connector ports, 46
 placement, 46

D

Delivery, 34
 box, 34
 Dimensions, 37
 instrument, 37
 Documentation, 8

E

Emergency, 16
 procedures, 16
 Environmental conditions, 39

F

FCC compliance, 115
 Fuse, 100
 replace, 100

I

Important user information, 5
 Inlet tubing, 66
 prime inlet tubing, 66
 prime inlets, 66
 Installation, 44, 54
 hardware, 44
 software, 54

Instrument control panel, 20, 21
 status indications, 21
 Instrument overview, 19, 20
 exterior design, 19
 instrument control panel, 20

L

Log on, 59
 UNICORN, 59

M

Maintenance, 93–95
 CIP, 95
 clean the instrument, 94
 Manufacturing information, 110

N

Notes and tips, 6

O

On site service, 121

P

Prepare sample for loading, 84, 85
 using a sample loop, 84
 using the pump, 85
 Prepare the system, 48, 59, 66
 prime inlets, 66
 purge system pumps, 66
 start UNICORN, 59
 waste tubing, 48
 Prime inlet tubing, 66
 Procedures after run, 90–92
 column CIP, 90
 column storage, 91
 shutting down, 92
 system storage, 91
 Procedures after the run, 90
 create CIP method, 90
 Product return or servicing, 122
 Pump, 66
 purge system pump, 66

Pump piston rinsing system, 50
 illustration, 50
 prime, 50
Pumps, 66
 purge system pumps, 66
Purpose of this manual, 6

R

Recycling information, 108
 decontamination, 108
 disposal of electrical components, 108
 recycling of hazardous substances, 108
Reference information, 105
 system specifications, 105
Regulatory information, 109
Run a method, 80, 81, 86–88
 create a method, 81
 evaluate the run, 88
 monitor/interact with the run, 87
 start a method run, 86

S

Safety notices, 11
Safety precautions, 11
 introduction, 11
Software, 28
 software modules, 28
Storage, 34
System Control, 29
 user interface, 29
System Control module, 30
 icons, 30
System recommendations, 42
 computer specifications, 42

T

Technical specifications, 105
 system specifications, 105
Troubleshooting, 102
Typographical conventions, 6

U

UK, 112
 conformity, 112

UKCA, 112
 marking, 112
UNICORN , 28, 59, 60
 connect to system, 60
 log on, 59
 software modules, 28
 start, 59

W

Waste tubing, 48
 prepare, 48

Page intentionally left blank

**Give feedback on this document**

Visit 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.

ÄKTA, ÄKTA avant, ÄKTA go, ÄKTA pcc, ÄKTA pure, Superloop, and UNICORN are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Decon and Decon 90 are a trademarks of Decon Laboratories Inc. Tween is a trademark of the Croda Group of Companies. Windows is a trademark of Microsoft group of companies.

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

© 2020–2023 Cytiva

UNICORN © 2020–2023 Cytiva

Any use of software may be subject to one or more end user license agreements, a copy of, or notice of which, are available on request.

For local office contact information, visit cytiva.com/contact

29360951 AE V:14 05/2023