

UNICORN™ software

Quick installation guide

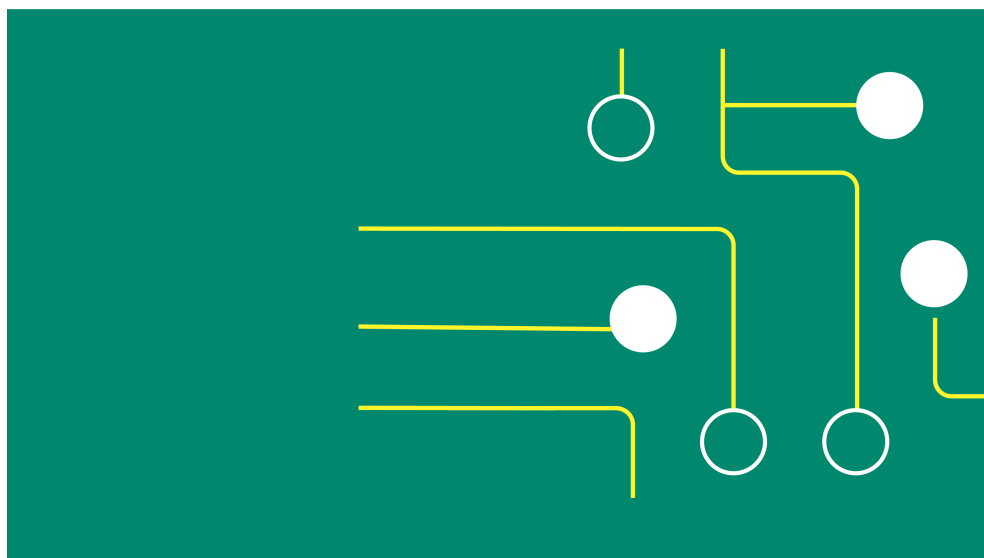


Table of Contents

1	Installation overview	3
2	Full installation	7
2.1	Step 1: Download the UNICORN software package	9
2.2	Step 2: Install the UNICORN software components	11
2.3	Step 3: Configure e-Licenses	17
2.4	Step 4: Configure the IP	23
2.5	Step 5: Define a system	31
3	Custom installation	34
3.1	Step 1: Download the UNICORN software package	36
3.2	Step 2: Install the UNICORN Database	38
3.3	Step 3: Install the Software Licensing Server	46
3.4	Step 4: Install the UNICORN software	55
3.5	Step 5: Configure the IP	62
3.6	Step 6: Define a system	70
4	Troubleshooting	71

1 Installation overview

About this document

This document contains an overview of the UNICORN™ software installation instructions. The screenshots included in this document might show slight differences with certain UNICORN versions and with certain Windows versions without affecting the instruction steps. For more comprehensive information about the UNICORN software installation, see the *UNICORN Administration and Technical Manual*.

About UNICORN

The UNICORN software can be installed on a computer to control an ÄKTA™ chromatography system connected to that computer. The software has the following components, which are installed one after another.

- UNICORN Database
- UNICORN Software
- Software Licensing Server

Note: *This document focuses on ÄKTA chromatography systems, but the UNICORN software can also be installed to control other types of systems such as for cultivation, filtration, and oligosynthesis.*

General computer specifications

For information on computer specifications, see the *UNICORN Computer specifications* document available on the Cytiva web.

For information about compatibility between UNICORN versions, the supported operating systems, database versions, and instrument configuration, see the UNICORN compatibility matrix at

<https://www.cytiva.com/UNICORNcompatibility>.

The following applies:

- A screen resolution of 1280x1024 or higher is recommended. Parts of the UNICORN user interface can not be displayed properly using a lower resolution.
- Changing the default font and changing the font size from 100% in Windows can cause problems in the UNICORN user interface.
- Windows power save features must be turned off to avoid conflicts with system operations.
- UNICORN is not compatible with the Windows feature High DPI Awareness, which allows the graphical user interface to be scaled. The interface scale must remain at 100% to avoid issues with clipping and misaligning of parts of the UNICORN user interface. Normally, the scale is set at 100% by default.

Installation types

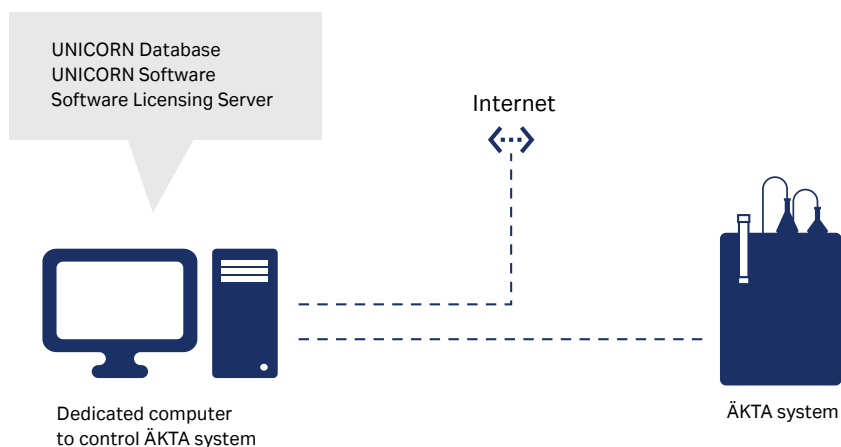
Depending on whether your computer is only connected to one ÄKTA system or connected to other computers controlling ÄKTA systems in a network, the UNICORN software can be installed in the following ways:

- Full installation
- Custom installation



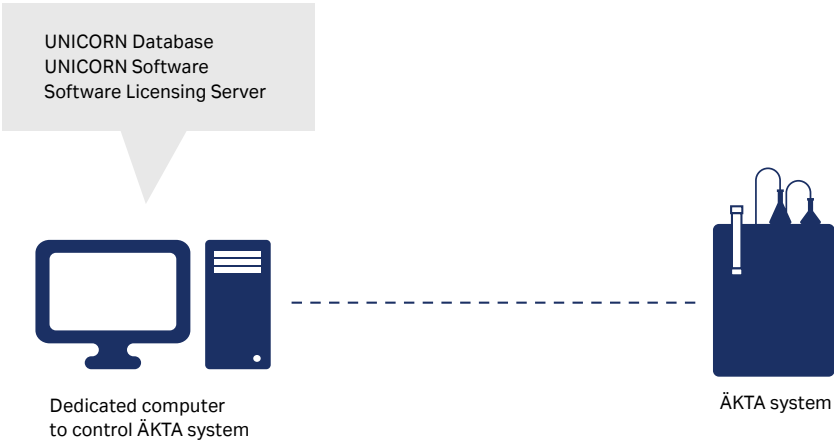
IMPORTANT

If you need to connect your dedicated computer to the Internet while connected to the ÄKTA system as in the illustration below, **you need two Network Interface Cards (NICs) installed in your dedicated computer.**



Full installation overview

When you have only one ÄKTA system, select the **Full installation** option. All the UNICORN software components are installed in the dedicated computer, see the illustration below.

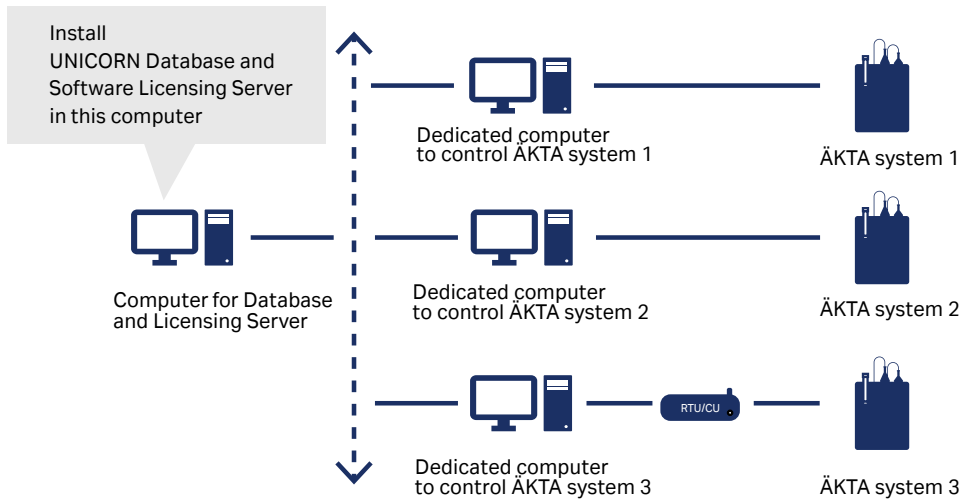


Custom installation overview

Select the **Custom installation** option when you have:

- **several ÄKTA systems,**
- all the dedicated computers connected to the same network, and
- an additional computer connected to the same network. This computer is where the **UNICORN software Database and Software Licensing Server** will be installed.

Note: You can choose to install the **UNICORN Database** on one computer and the **License Server** on another.



2 Full installation



IMPORTANT

You need an **Activation ID** to download the UNICORN software installation package and to retrieve an e-License from the Cytiva software e-Licensing web site. The activation ID together with the information about the software is sent via e-mail. Contact your Cytiva representative if you have not received an activation ID.

An Internet connection is needed to download the UNICORN software installation package and to retrieve the e-License.

UNICORN Database
UNICORN Software
Software Licensing Server



Dedicated computer
to control ÄKTA system



ÄKTA system

In this chapter

Section		See page
2.1	Step 1: Download the UNICORN software package	9
2.2	Step 2: Install the UNICORN software components	11
2.3	Step 3: Configure e-Licenses	17
2.4	Step 4: Configure the IP	23
2.5	Step 5: Define a system	31

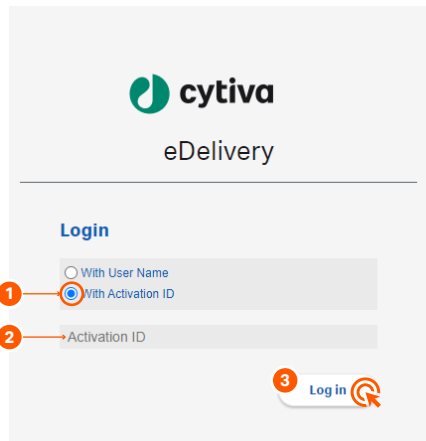
2.1 Step 1: Download the UNICORN software package



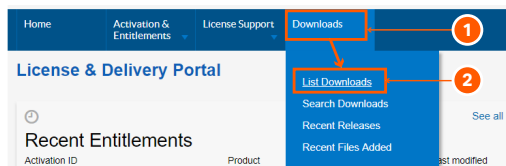
IMPORTANT

Make sure that all the Windows updates are completed. Otherwise, the computer might start installing Windows updates during the UNICORN software installation and significantly delay the installation process.

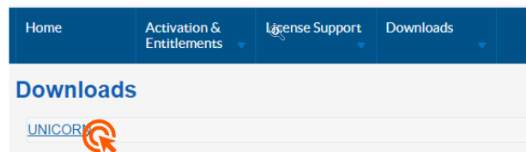
Step	Action
1	Enter the following address: https://www.cytiva.com/eDelivery .
2	Follow the steps in the illustration below to log in.



3 Follow the steps in the illustration below.

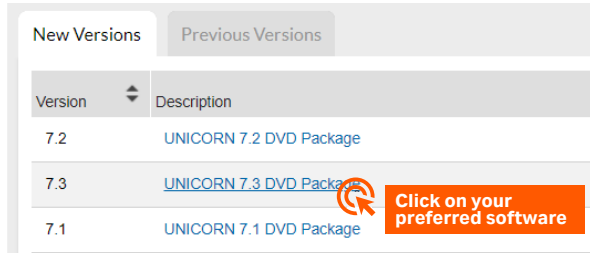


4 Enter the UNICORN page.

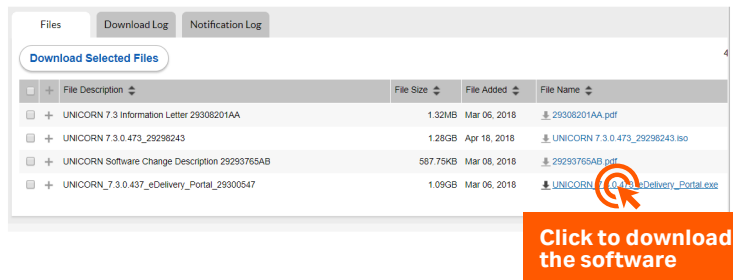


Step	Action
------	--------

5	Select a software version.
---	----------------------------



6	Download the software.
---	------------------------



The software package is downloaded in the **Downloads** folder.

7	Go to Section 2.2 Step 2: Install the UNICORN software components, on page 11.
---	--

2.2 Step 2: Install the UNICORN software components

It is assumed that you have downloaded the UNICORN software installation package on the computer where the UNICORN software will be installed.

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

1	On the computer where the UNICORN software installation package is downloaded, locate the downloaded installation package.
---	--

2	Double-click the UNICORN_7.x.x.xxx_eDelivery_Portal.exe file.
---	--

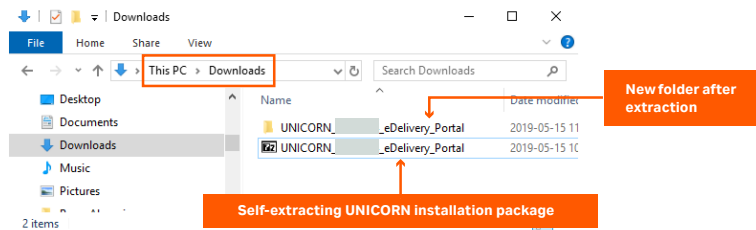


IMPORTANT

The extraction can take **several minutes**.

If the **Open File - Security Warning** window appears, click **Run**.

3	The self-extracting function creates a folder with the same name as the downloaded file at the same location.
---	---



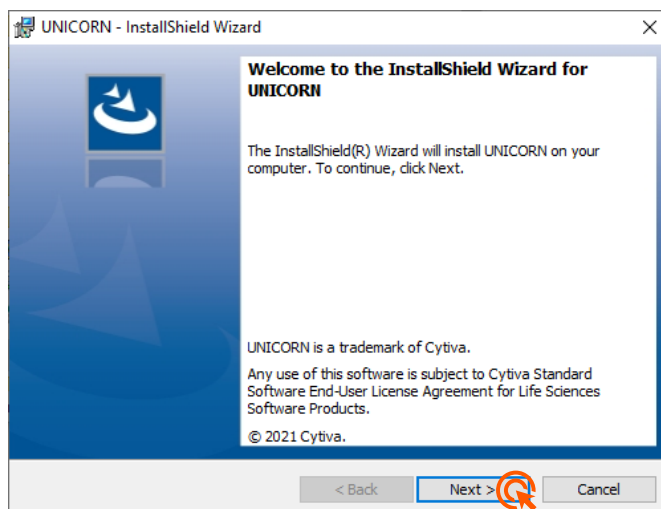
Double-click the **UNICORN_7.x.x.xxx_eDelivery_Portal** folder.

4	Double-click Setup.exe .
---	---------------------------------

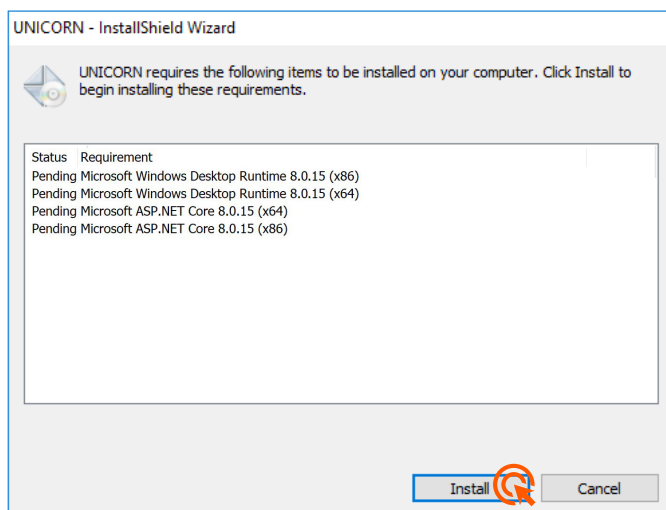
If the **Open File - Security Warning** window appears, click **Run**.

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

5	Click Next .
---	---------------------

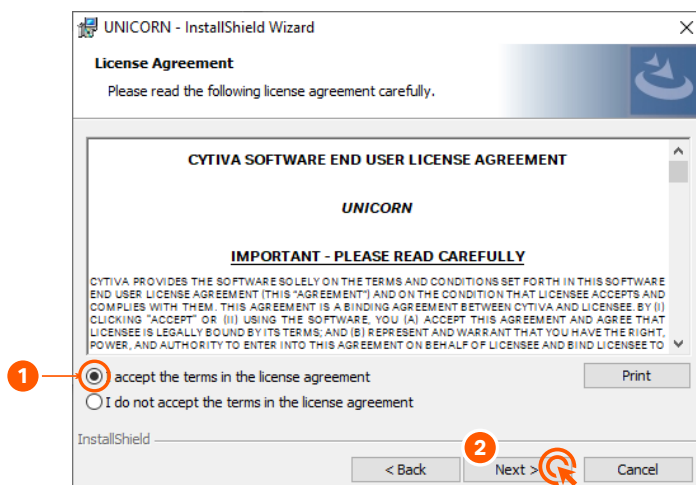


6	If your computer is missing any software requirements , the installer will install them. The computer can restart several times. After the restart, double-click the Setup.exe again, see step 5 .
---	--

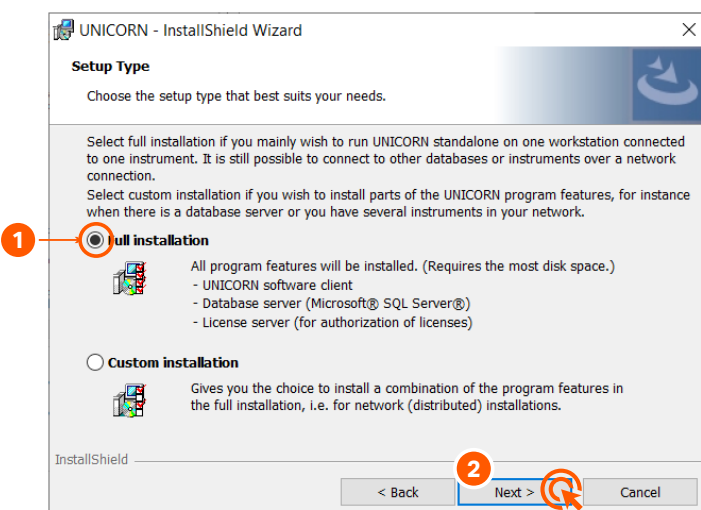


Step	Action
------	--------

7	Accept the license agreement and click Next .
---	--

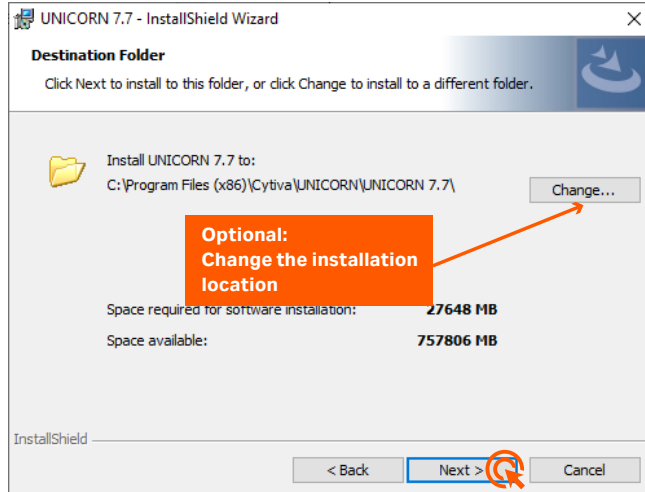


8	Select the installation type and click Next .
---	--

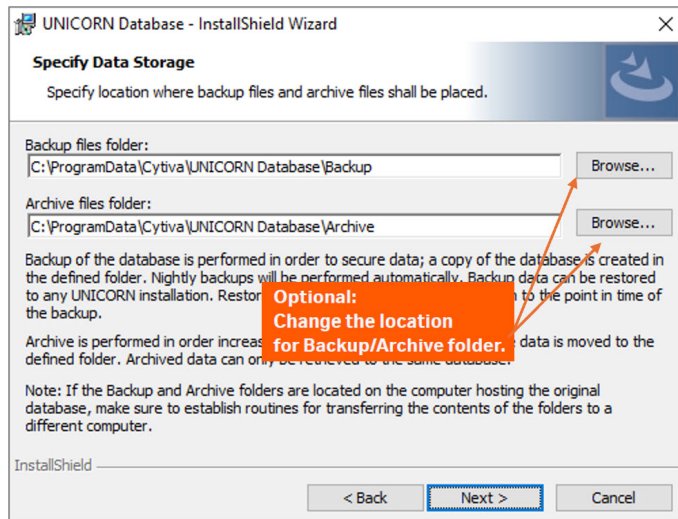


Step	Action
------	--------

9	Click Next .
---	---------------------

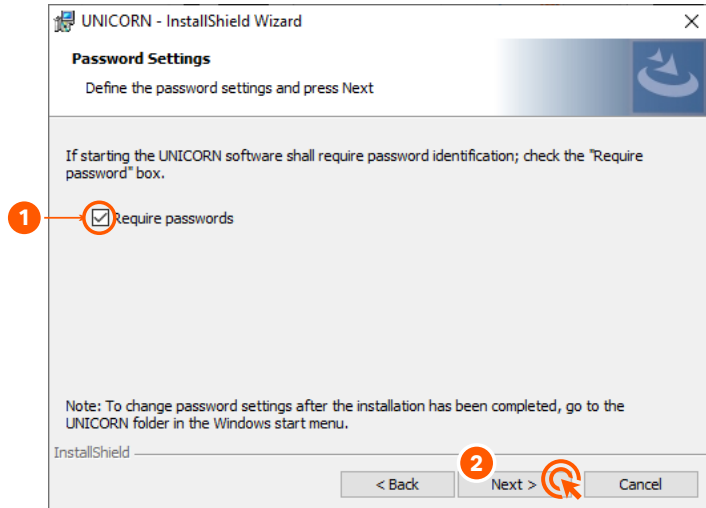


10	Click Next .
----	---------------------



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

11	Select Require password (optional).
----	--



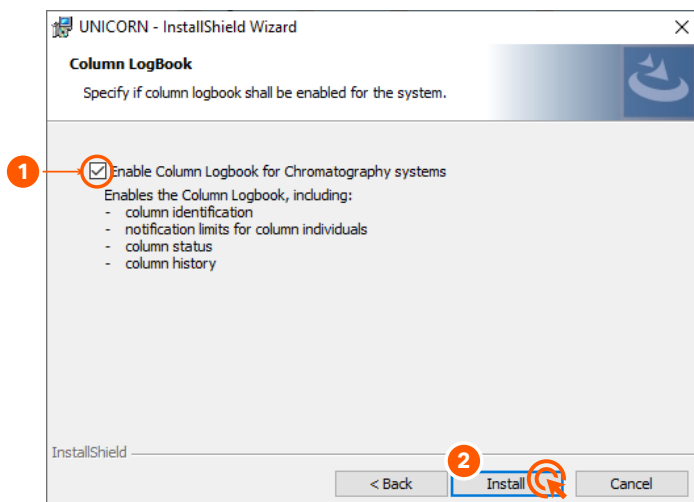
12	
----	--

**IMPORTANT**

Enable **Column Logbook** only if you have a **license** for it.

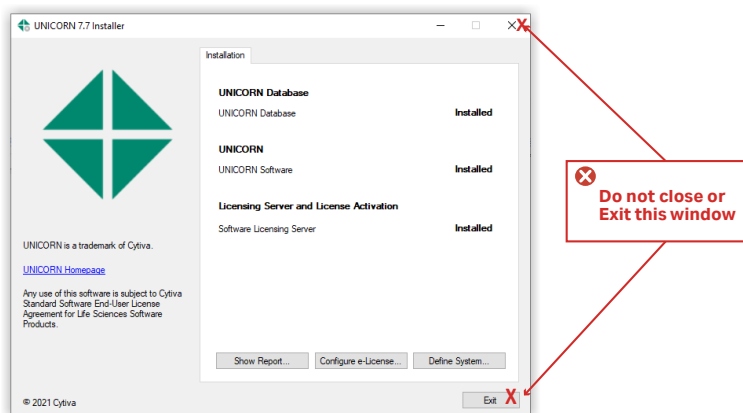
Step	Action
------	--------

	Select Column Logbook and then click Install .
--	--



Wait, this step can take several minutes.

If the following dialog appears, the first step for the installation is complete.



IMPORTANT

Do not close the UNICORN Installer window (see the illustration above). If you accidentally close it, perform step 4 of this procedure again.

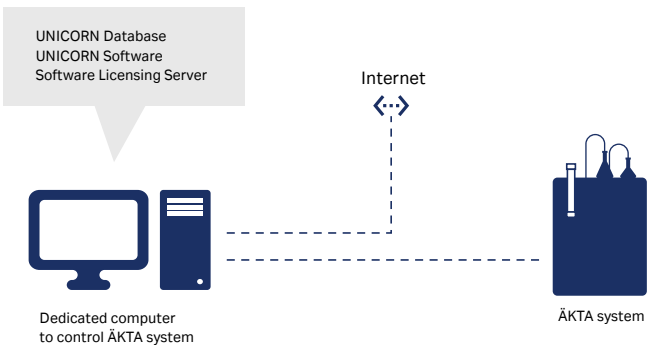
2.3 Step 3: Configure e-Licenses



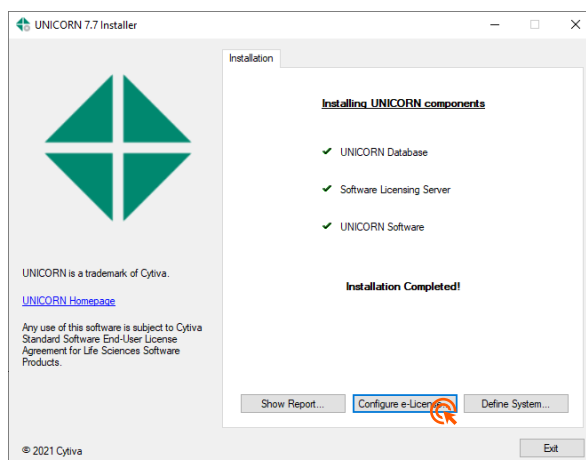
IMPORTANT

To configure an e-License, **you must have the Activation ID** and an Internet connection.

If the dedicated computer is connected to the Internet and the ÄKTA system as illustrated below, you can retrieve the e-License using the dedicated computer. Make sure that the dedicated computer has an Internet connection.

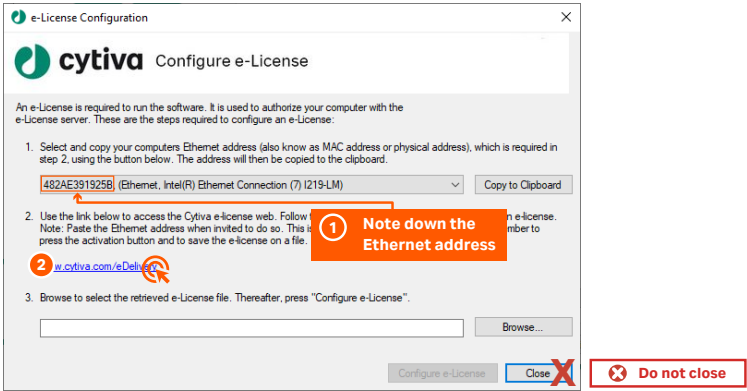


Step	Action
1	Open the Configuration e-License window as shown below.



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

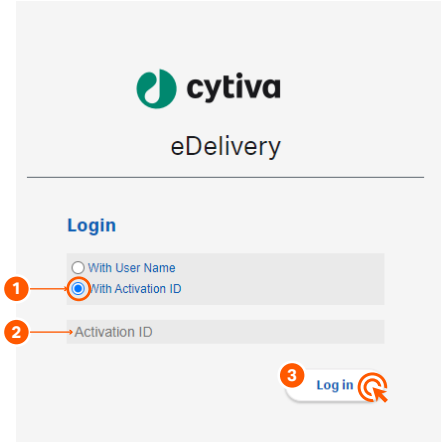
2	Follow the steps in the illustration below to configure an e-License.
---	---



IMPORTANT

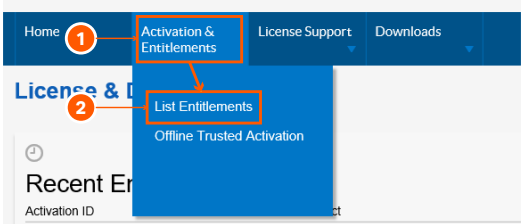
Note down your Ethernet address. The Ethernet address stated in the illustration above is just **an example**.

3	Follow the steps in the illustration below to log in.
---	---

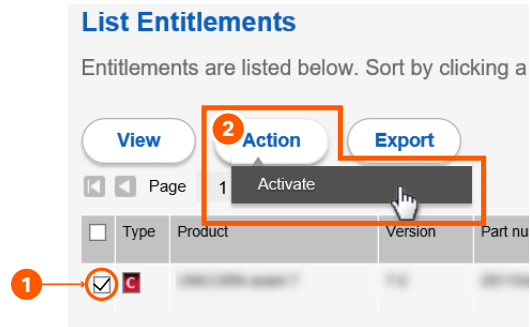


Step Action

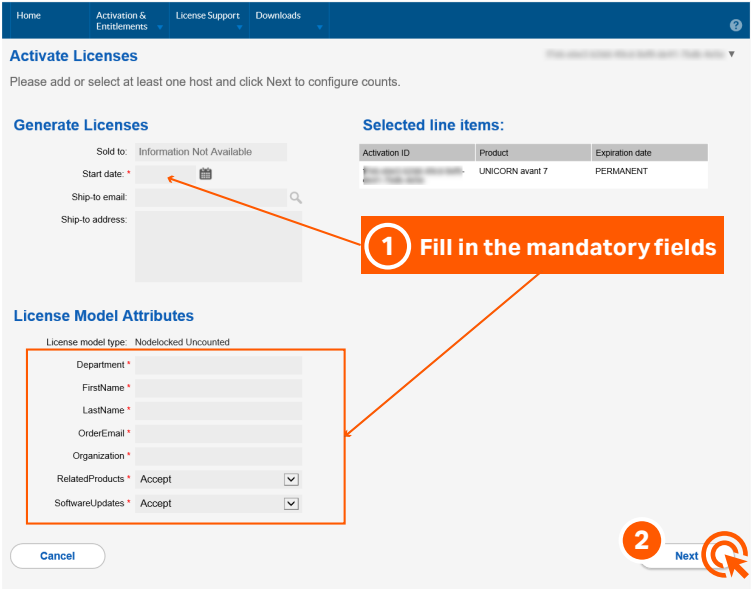
4 Follow the steps in the illustration below.



5 Follow the steps in the illustration below.



6 Follow the steps in the illustration below.



Step Action

7 Follow the steps in the illustration below.

Configure Hosts

Please enter values for the hosts and click Verify to validate the license. Or click Generate to generate the license.

Configure Hosts

Server hosts: (no license hosts selected)

Selected line items:

Activation ID	Product	Expiration date
UNICORN 7.0	UNICORN avant 7	PERMANENT

Create License Host

Please select a license host type and enter an appropriate value.

Node-locked hosts: ETHERNET =

2 Enter the Ethernet address you have noted in step 2.

Configure Hosts

Please enter values for the hosts and click Verify to validate the license. Or click Generate to generate the license.

Configure Hosts

Server hosts: A4C3F0B51775

Selected line items:

Activation ID	Product	Expiration date
UNICORN 7.0	UNICORN avant 7	PERMANENT

License details:

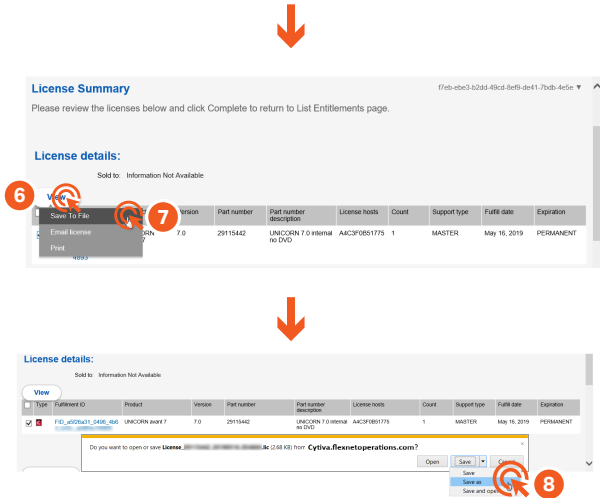
Sold to: Information Not Available

View

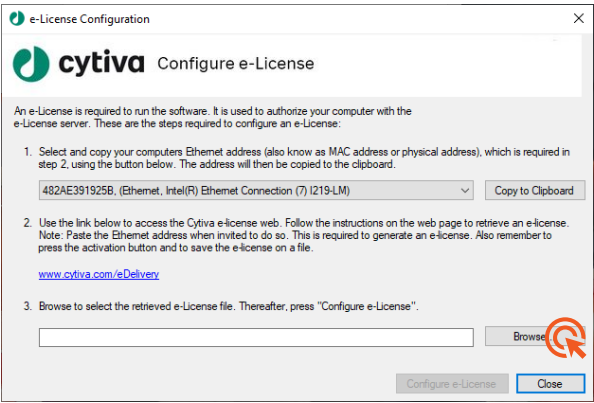
Type	Fulfillment ID	Product	Version	Part number	Part number description	License hosts	Co
☒	FID_a5f26a31_04f6_4b64_b43c_a5964e1f4893	UNICORN avant 7	7.0	29115442	UNICORN 7.0 internal no DVD	A4C3F0B51775	1

Step Action

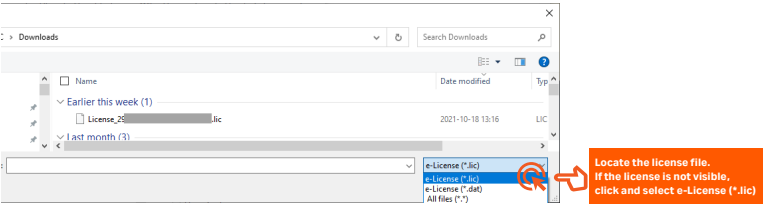
8 Follow the steps in the illustration below.



9 Go back to the **e-License Configuration** window and click **Browse**.



10 Browse to the e-License location.



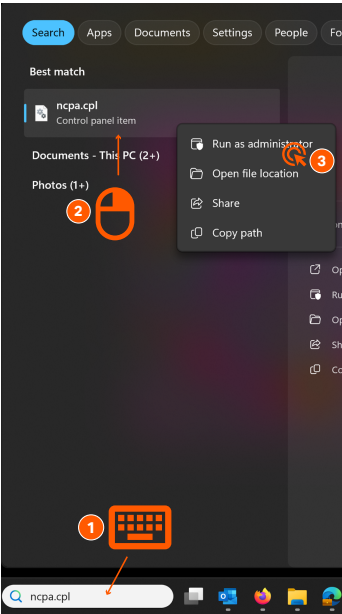
2.4 Step 4: Configure the IP

Step	Action
1	Make sure that your dedicated computer is connected to the Internet and to the ÄKTA system.
2	Click the Windows search icon.



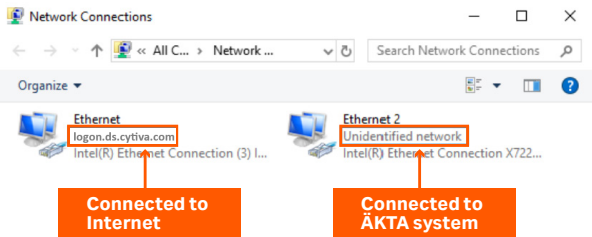
Step Action

3 Type ***ncpa.cpl*** in the **search box**, and then press **Enter**.



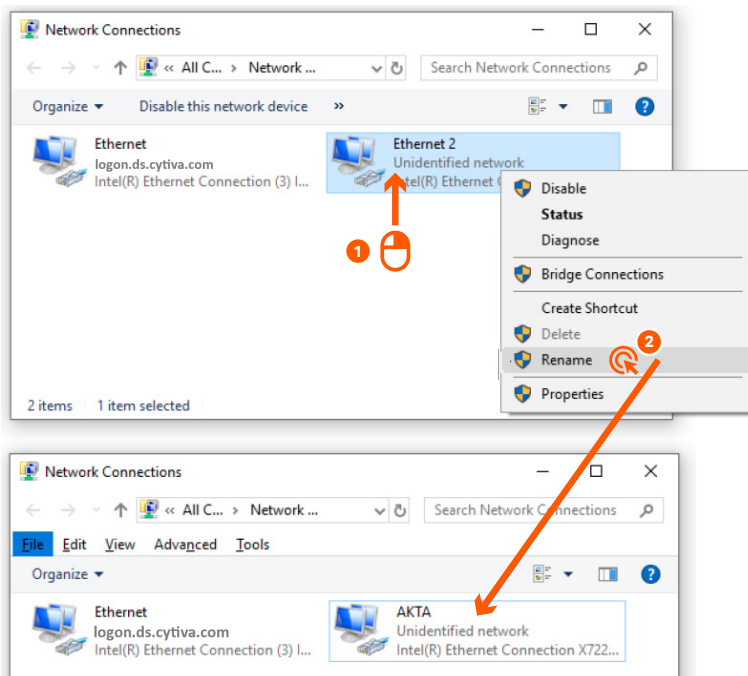
IMPORTANT

If you have two Ethernet cards, one Ethernet card is connected to the Internet (**Ethernet in the illustration below**) and the other (**Ethernet 2 in the illustration below**) is connected to the ÄKTA system. Generally, the Ethernet card connected to the Internet has the company domain name. For example, in the illustration below, it is **logon.ds.ge.com**. The Ethernet card connected to the ÄKTA system generally states **Unidentified network** as seen in the illustration below.

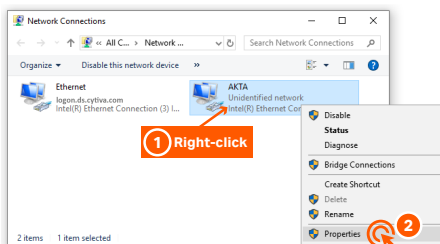


Step	Action
------	--------

- | | |
|---|--|
| 4 | Rename the Ethernet 2 connection to AKTA . |
|---|--|

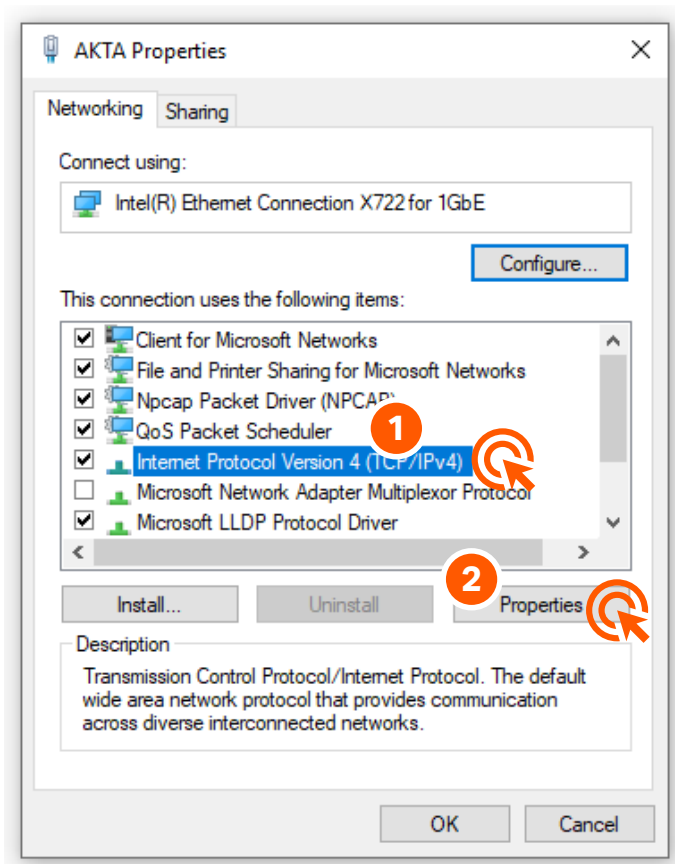


- | | |
|---|---|
| 5 | Right-click the AKTA network connection and select Properties . |
|---|---|



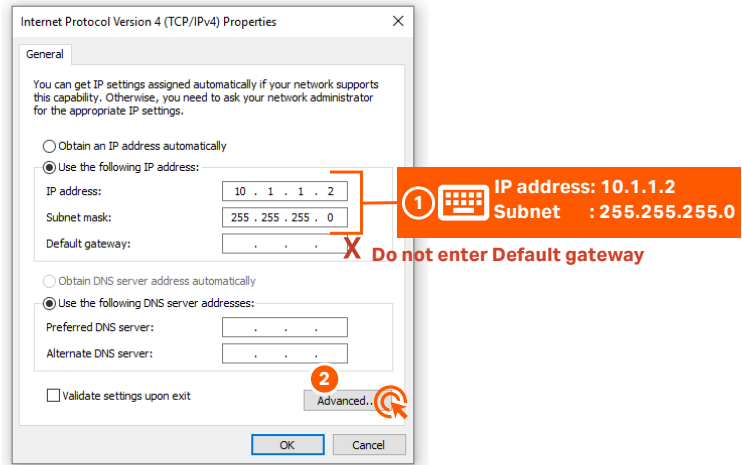
Step	Action
------	--------

6	Select Internet Protocol Version 4 .
---	---

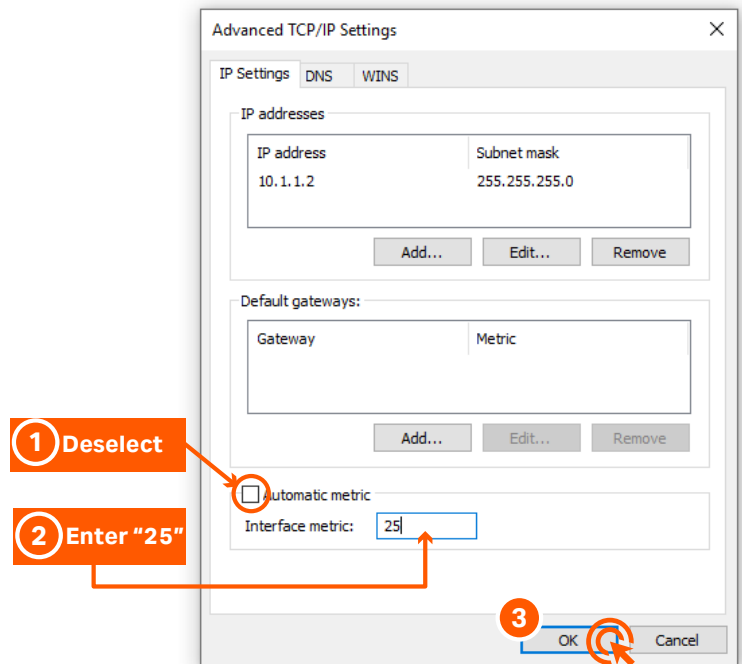


Step	Action
------	--------

- | | |
|---|--|
| 7 | Enter the IP address and the Subnet mask as shown in the illustration below. |
|---|--|



- | | |
|---|--|
| 8 | Apply the advanced TCP/IP settings as shown in the illustration below. |
|---|--|



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

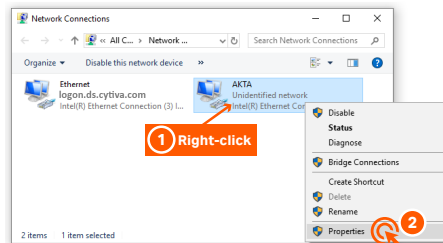
9	Click OK or Close in all the open dialogs.
---	--



IMPORTANT

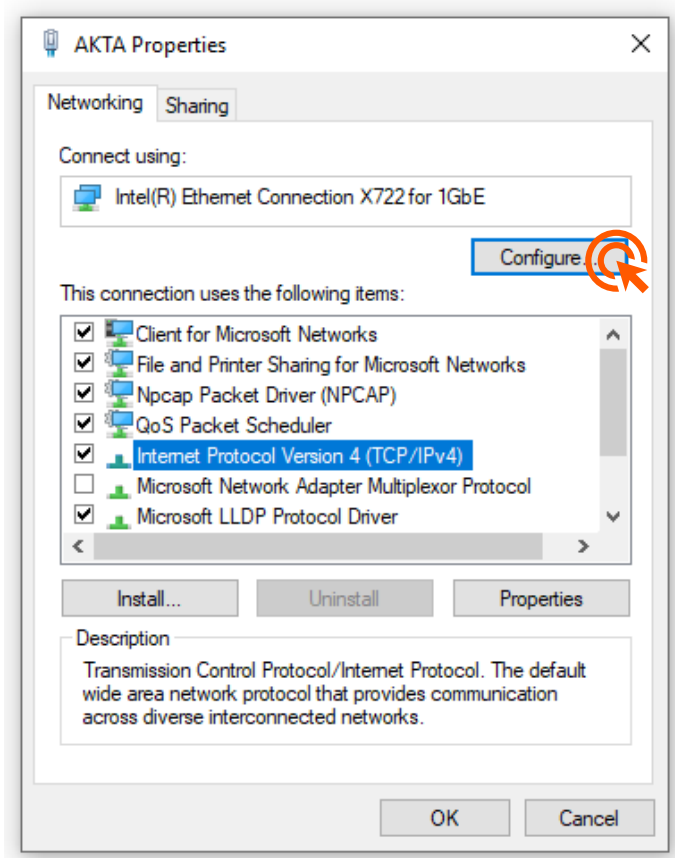
- If the system to be connected is a CU-based system, for example, ÄKTAprouess™, ÄKTA ready™, ÄKTA readyflux™, etc., perform steps 10 to 12.
- If the system to be connected is not a CU-based system, for example, ÄKTA avant, ÄKTA pure™, ÄKTA go™, ÄKTA pilot™ 600, ÄKTA process™, etc., skip steps 10 to 12. The **Speed & Duplex** property for these systems is set to **Auto Negotiation** by default.

10	Access the Properties window of the ÄKTA network connection.
----	---



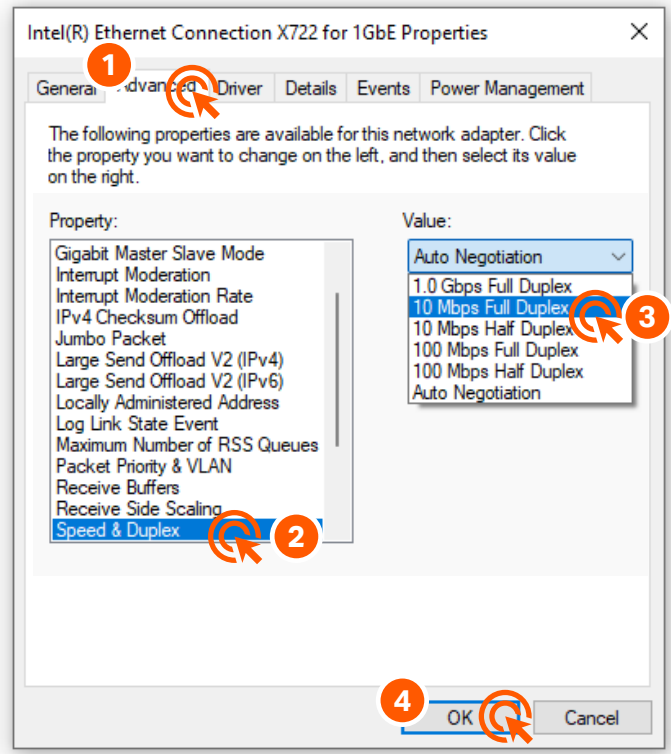
Step	Action
------	--------

11	Click Configure...
----	---------------------------



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

- | | |
|----|--|
| 12 | Set the Speed & Duplex property value to 10 Mbps Full Duplex as shown in the illustration below. |
|----|--|



- | | |
|----|---|
| 13 | Go to Section 2.5 Step 5: Define a system, on page 31 . |
|----|---|

2.5 Step 5: Define a system



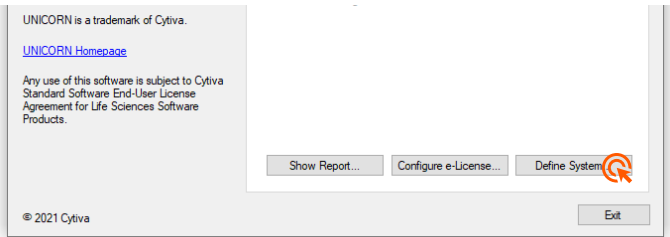
IMPORTANT

You need an **Instrument Configuration (IC)** to define your system.

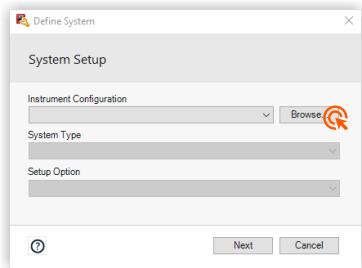
The latest IC can be downloaded from the product page of your system. Go to <https://www.cytiva.com/akta> and search for your ÄKTA system on the page. The IC can be found under **RELATED DOCUMENTS → SOFTWARE** on your ÄKTA system web page.

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

- | | |
|---|---------------------------------------|
| 1 | Open the Define System window. |
|---|---------------------------------------|



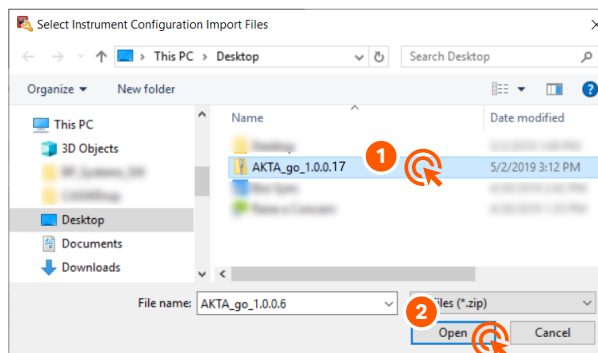
- | | |
|---|--|
| 2 | Browse to the Instrument Configuration location. |
|---|--|



Click and browse to the location where the Instrument Configuration is downloaded and unpacked.

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

- | | |
|---|---|
| 3 | Open the Instrument Configuration file. |
|---|---|

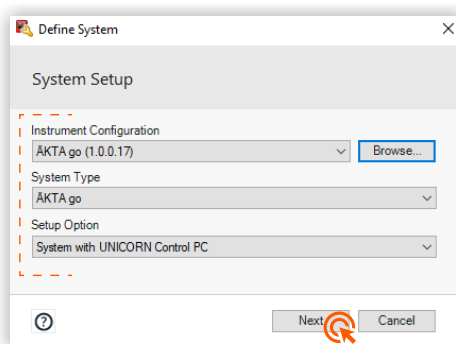


Note:

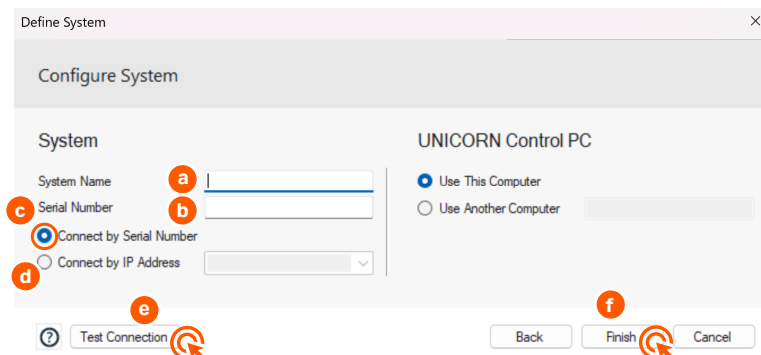
Both **.zip** and **.ic** are valid Instrument Configuration extensions.

- | | |
|---|----------------------------------|
| 4 | Select the system setup options. |
|---|----------------------------------|

These fields are automatically populated with default values. Use the drop-down list to change.



- | | |
|---|---|
| 5 | Follow the steps below to configure the system. |
|---|---|



Step	Action
	a. Enter your preferred system name. This name cannot be changed later. b. Enter your ÄKTA serial number. The number can be found on a sticker at the back of the system. c. Select Connect by Serial Number. d. Alternatively, you can select Connect by IP Address and enter the corresponding IP address. e. Click Test Connection. If the connection is successful, proceed to the next step. f. Click Finish.
6	When you have successfully defined your system, you can close the Define System dialog. Now the ÄKTA system is ready for use.

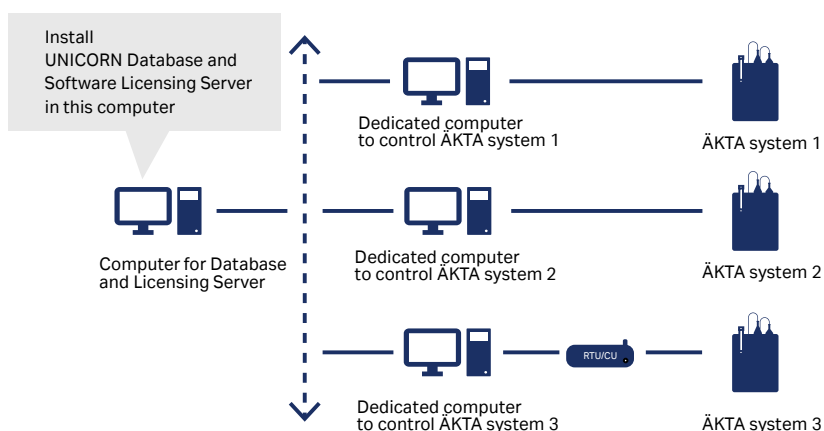
3 Custom installation



IMPORTANT

You need an **Activation ID** to download the UNICORN software installation package and to retrieve an e-License from the Cytiva software e-Licensing web site. The activation ID together with the information about the software is sent via e-mail. Contact your Cytiva representative if you have not received an activation ID.

An Internet connection is needed to download the UNICORN software installation package and to retrieve the e-License.



In this chapter

Section		See page
3.1	Step 1: Download the UNICORN software package	36
3.2	Step 2: Install the UNICORN Database	38
3.3	Step 3: Install the Software Licensing Server	46
3.4	Step 4: Install the UNICORN software	55
3.5	Step 5: Configure the IP	62
3.6	Step 6: Define a system	70

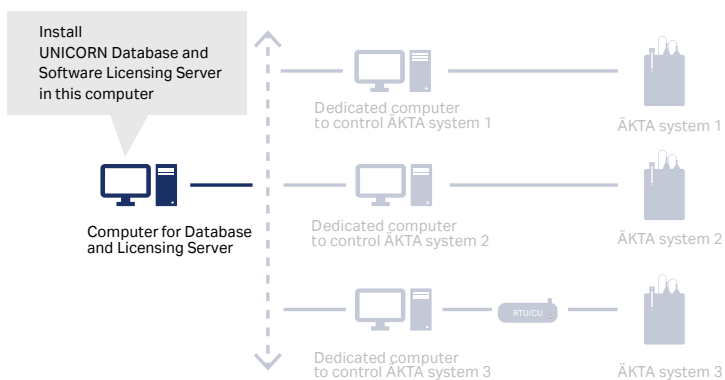
3.1 Step 1: Download the UNICORN software package



IMPORTANT

Make sure that all the Windows updates are completed. Otherwise, the computer might start installing Windows updates during the UNICORN software installation and significantly delay the installation process.

Step	Action
1	On the computer where the UNICORN software Database and the Software Licensing server will be installed, open any web browser.

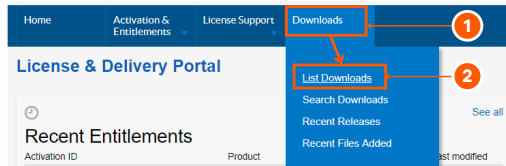


2 Enter the following address: <https://www.cytiva.com/eDelivery>.

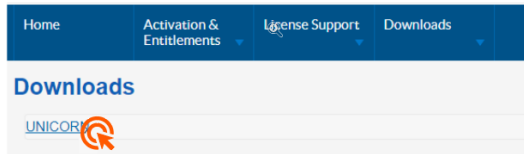
3 Follow the steps in the illustration below to log in.

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

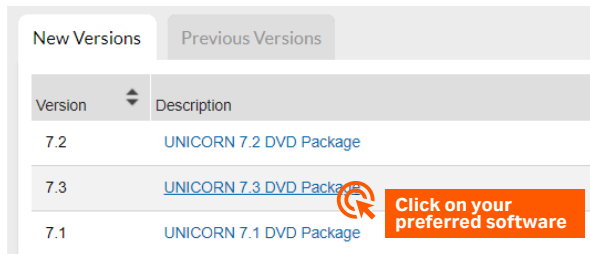
4	Follow the steps in the illustration below.
---	---



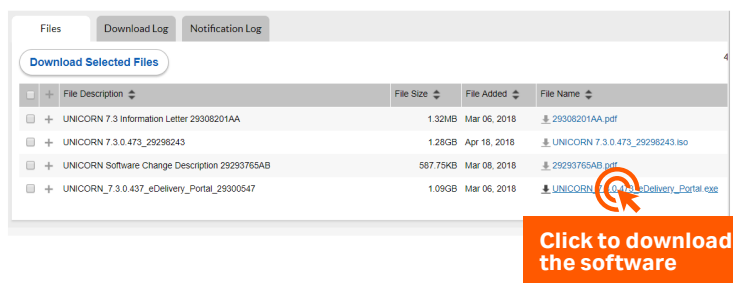
5	Enter the UNICORN page.
---	-------------------------



6	Select a software version.
---	----------------------------



7	Download the software.
---	------------------------



The software package will be downloaded in the **Downloads** folder.

8	Go to Section 3.2 Step 2: Install the UNICORN Database, on page 38 .
---	--

3.2 Step 2: Install the UNICORN Database

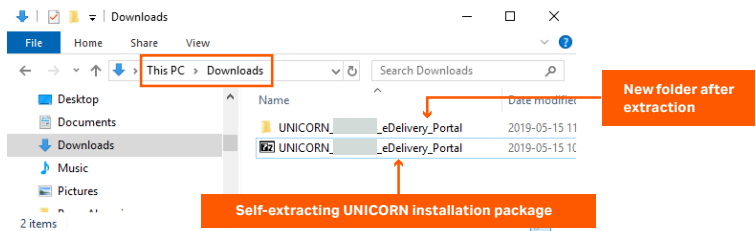
It is assumed that you have downloaded the UNICORN software installation package on the computer where the UNICORN software will be installed.

Step	Action
1	On the computer where the UNICORN software installation package is downloaded, locate the downloaded installation package.
2	Double-click the UNICORN_7.x.x.xxx_eDelivery_Portal.exe file.



IMPORTANT
The extraction can take **several minutes**.
If the **Open File - Security Warning** window appears, click **Run**.

- 3 The self-extracting function creates a folder with the same name as the downloaded file at the same location.



Double-click the **UNICORN_7.x.x.xxx_eDelivery_Portal** folder.

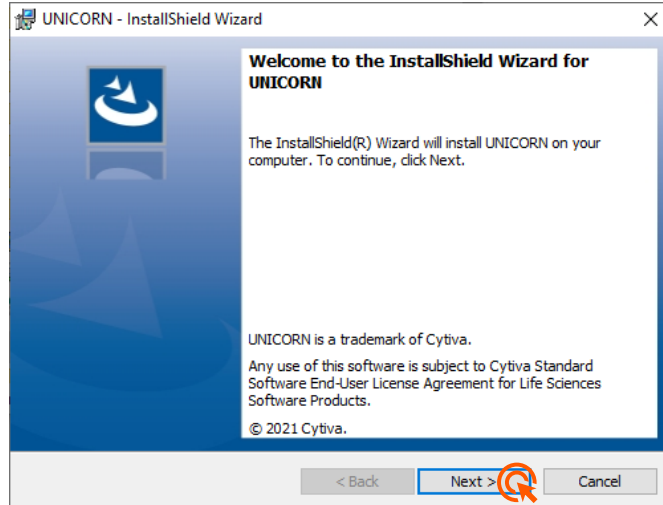


IMPORTANT
If your organization allows you to use **USB memory sticks**, copy the extracted installation package into a **USB memory stick** to simplify the installation process. **Otherwise**, make sure that the extracted files are accessible from all the related computers for this installation.

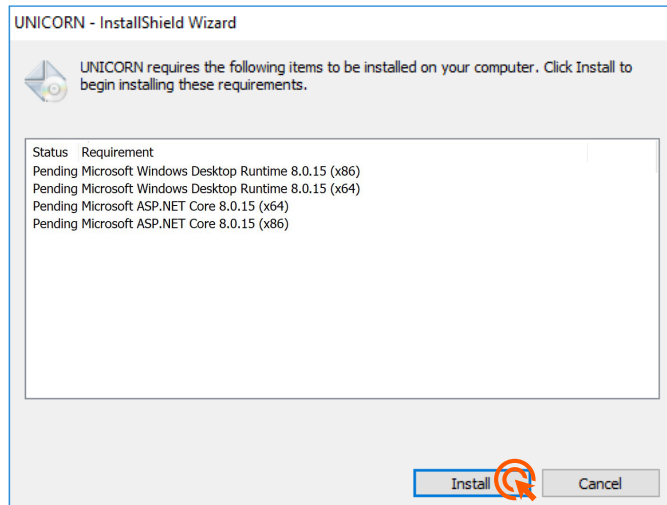
- 4 Double-click **Setup.exe**. If the **Open File - Security Warning** window appears, click **Run**.

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

- | | |
|---|---------------------|
| 5 | Click Next . |
|---|---------------------|

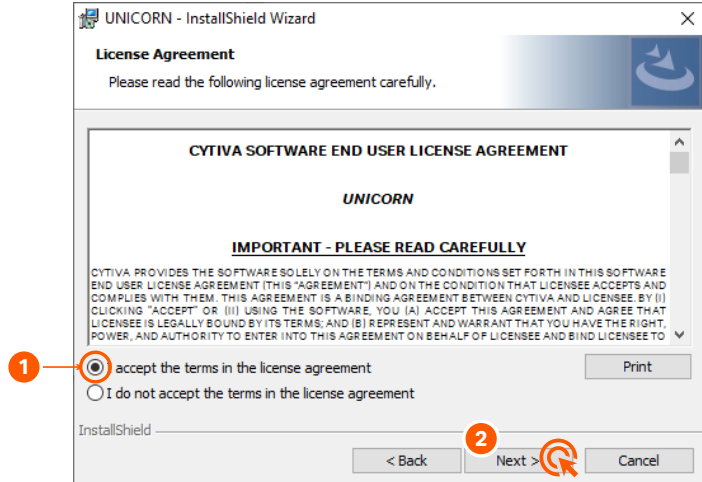


- | | |
|---|--|
| 6 | If your computer is missing any software requirements , the installer will install them. The computer can restart several times. After the restart, double-click the Setup.exe again, see step 5 . |
|---|--|

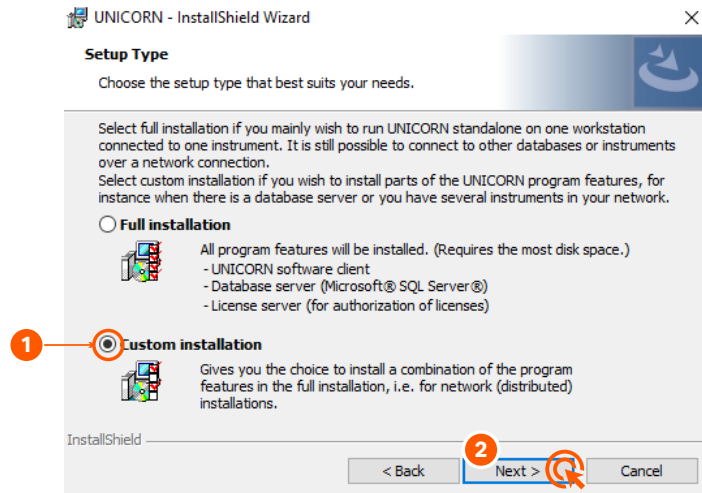


Step	Action
------	--------

7	Accept the license agreement and click Next .
---	--

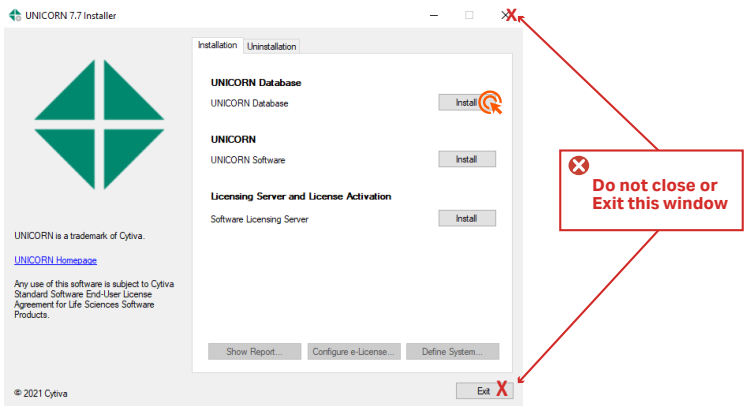


8	Select the installation type and click Next .
---	--

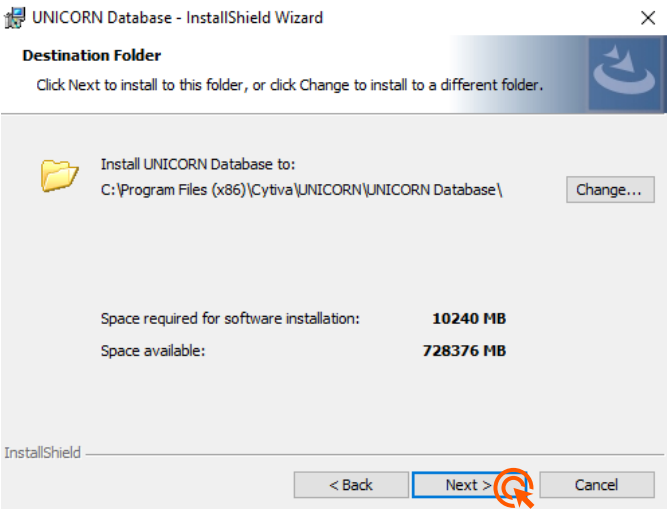


Step	Action
------	--------

9	Click Install next to UNICORN Database .
---	--

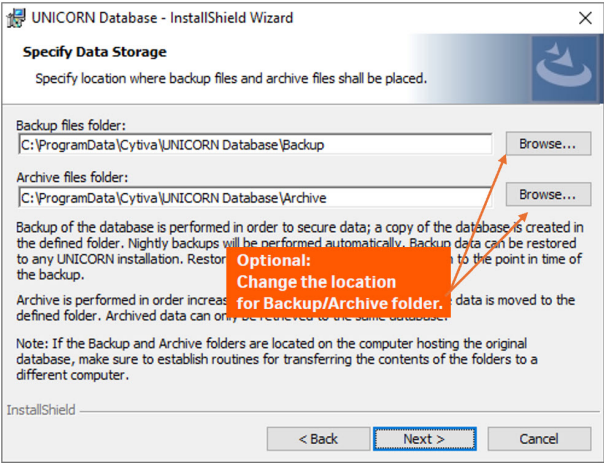


10	Click Next .
----	---------------------

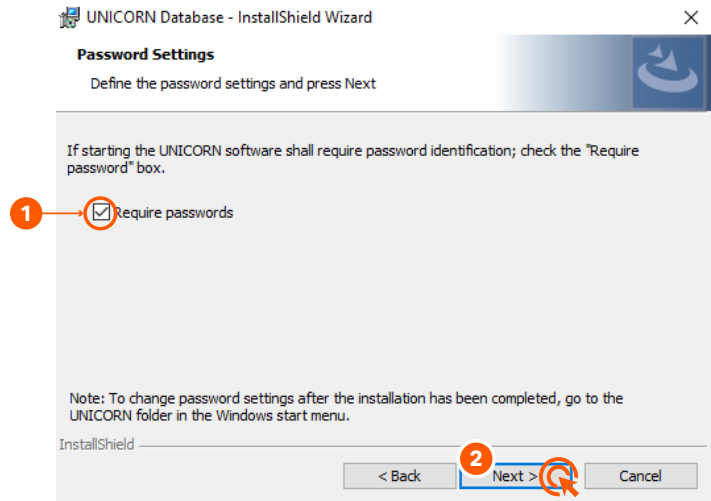


Step Action

11 Click **Next**.



12 Select **Require password** (optional).



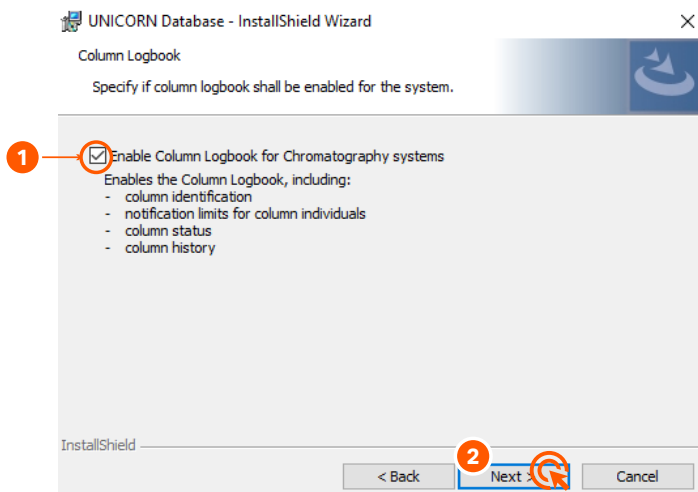
13

IMPORTANT

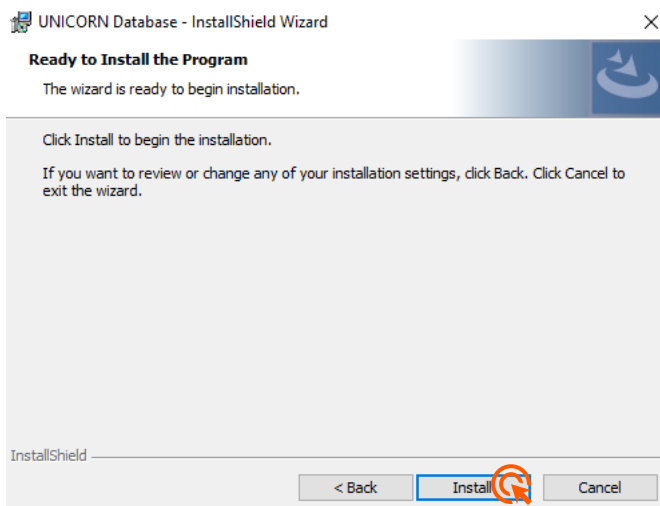
Enable **Column Logbook** only if you have a **license** for it.

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

Follow the steps in the illustration below.

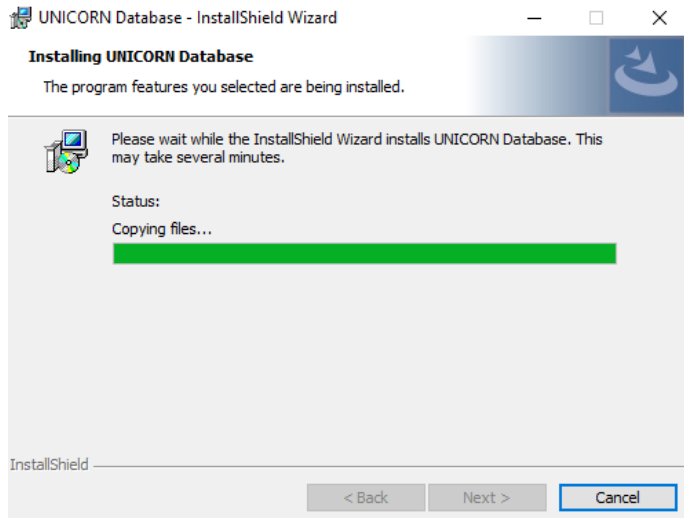


14 Click **Install**.

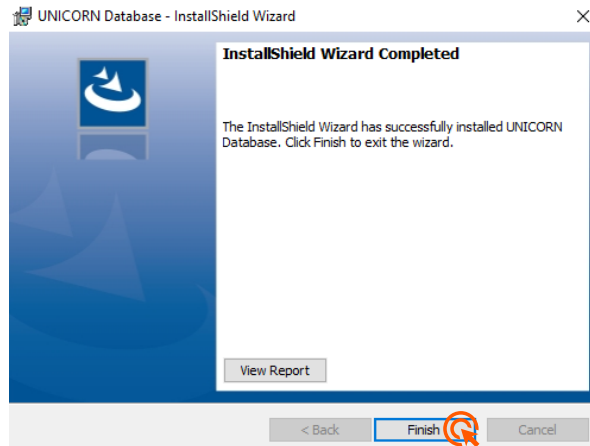


Wait.

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

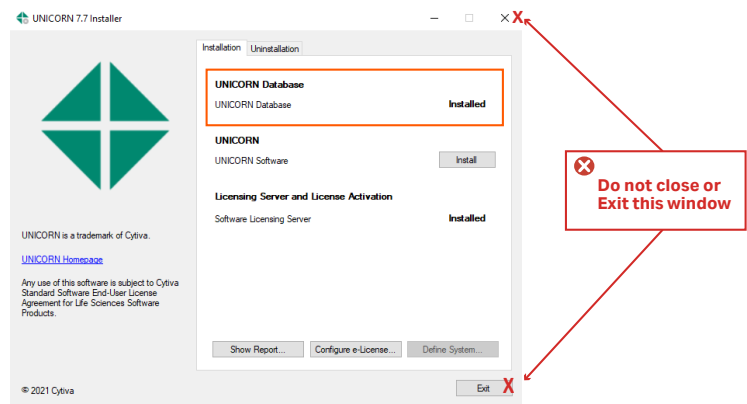


15 Click **Finish**.



If the following window appears, the database installation is finished.

Step Action



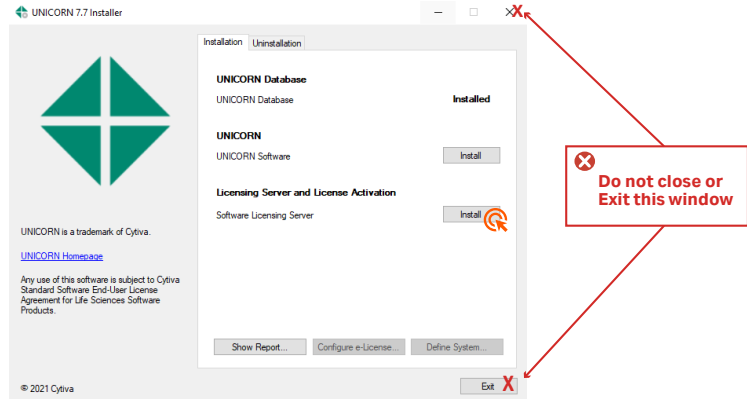
IMPORTANT

Do not close the UNICORN Installer window (see the illustration above). If you accidentally close it, perform steps 4 to 6 of this procedure again.

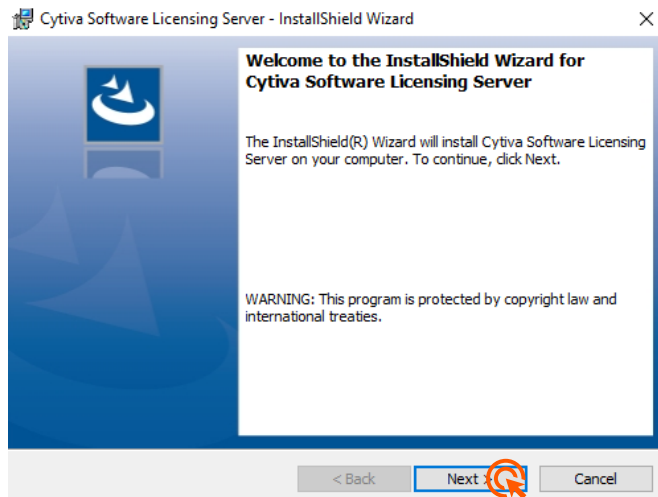
3.3 Step 3: Install the Software Licensing Server

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

- | | |
|---|---|
| 1 | Click Install next to Software Licensing Server . |
|---|---|

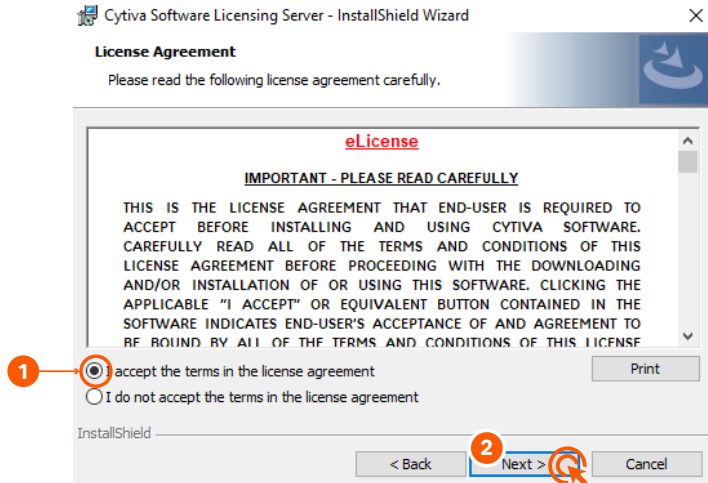


- | | |
|---|---------------------|
| 2 | Click Next . |
|---|---------------------|

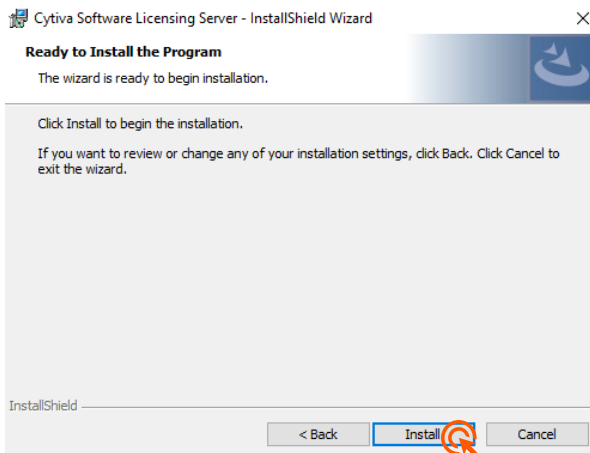


Step	Action
------	--------

3	Accept the license agreement and click Next .
---	--

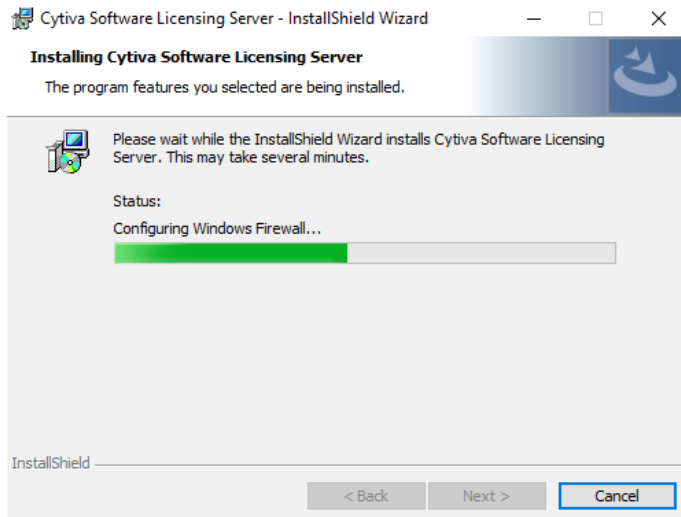


4	Click Install .
---	------------------------

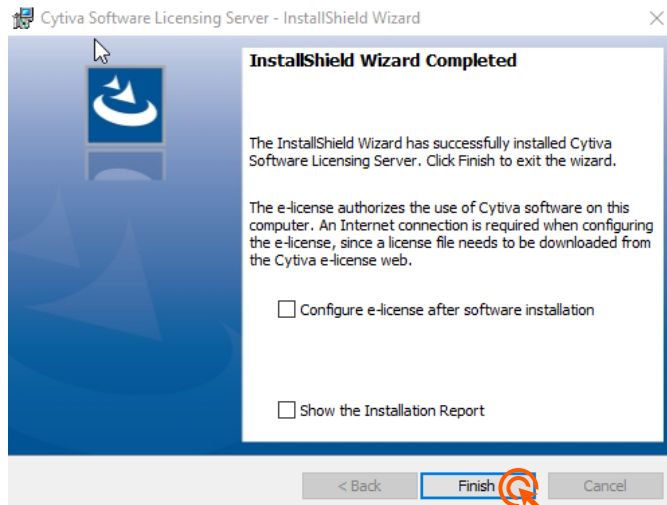


Wait.

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

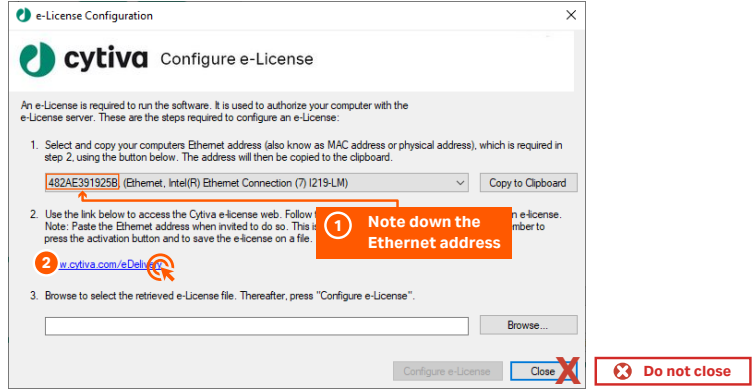


5 Click **Finish**.



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

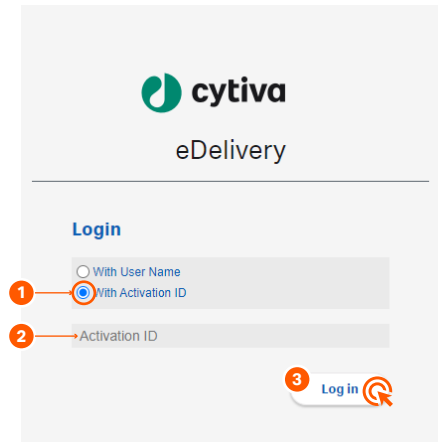
6	Follow the steps in the illustration below to configure an e-License.
---	---



IMPORTANT

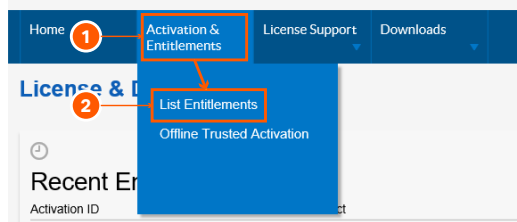
Note down your Ethernet address. The Ethernet address stated in the illustration above is just **an example**.

7	Follow the steps in the illustration below to log in.
---	---

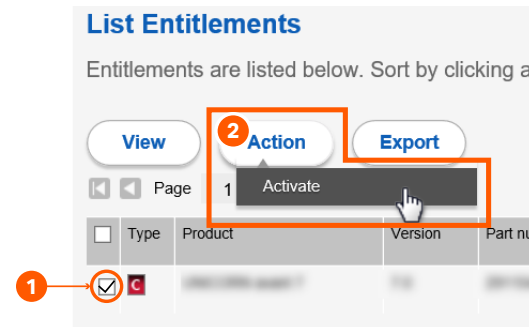


Step	Action
------	--------

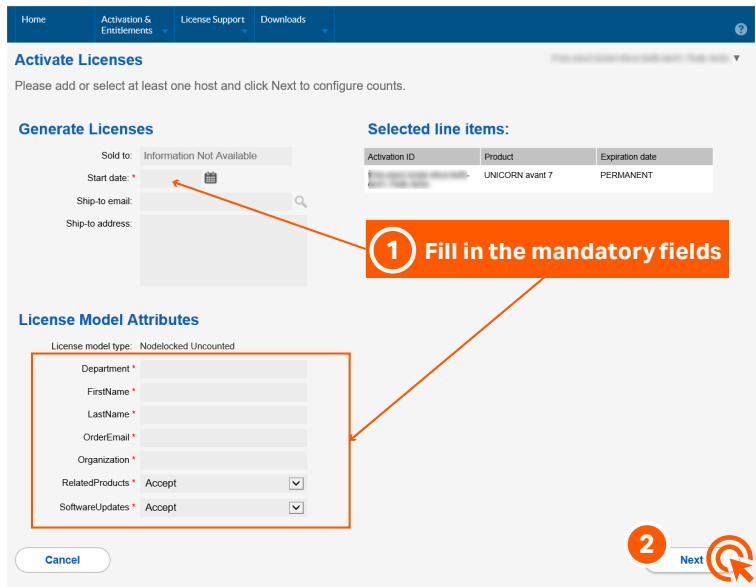
- | | |
|---|---|
| 8 | Follow the steps in the illustration below. |
|---|---|



- | | |
|---|---|
| 9 | Follow the steps in the illustration below. |
|---|---|



- | | |
|----|---|
| 10 | Follow the steps in the illustration below. |
|----|---|



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

11	Follow the steps in the illustration below.
----	---

Configure Hosts

Please enter values for the hosts and click Verify to validate the license. Or click Generate to generate the license.

Configure Hosts

Server hosts: (no license hosts selected)

Selected line items:

Activation ID	Product	Expiration date
...	UNICORN avant 7	PERMANENT

Buttons: Cancel, Back, Generate

Create License Host

Please select a license host type and enter an appropriate value.

Node-locked hosts: ETHERNET

Buttons: OK, Cancel

Callout: Enter the Ethernet address you have noted in step 2.

Configure Hosts

Please enter values for the hosts and click Verify to validate the license. Or click Generate to generate the license.

Configure Hosts

Server hosts: A4C3F0B51775

Selected line items:

Activation ID	Product	Expiration date
...	UNICORN avant 7	PERMANENT

Buttons: Cancel, Back, Generate

License details:

Sold to: Information Not Available

View

Type	Fulfillment ID	Product	Version	Part number	Part number description	License hosts	Co
☒	FID_a5f26a31_04f6_4b64_b43c_a5964e1f4893	UNICORN avant 7	7.0	29115442	UNICORN 7.0 internal no DVD	A4C3F0B51775	1

12 Follow the steps in the illustration below.



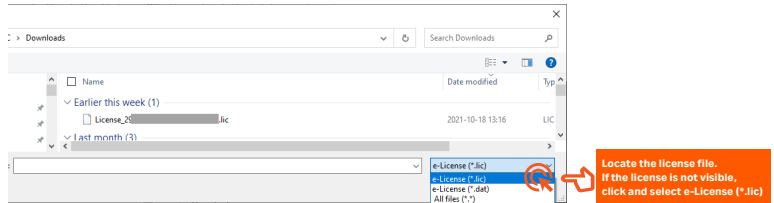
We recommend renaming the license so that the license can be easily identified later.

13 Go back to the **e-License Configuration** window and click **Browse**.

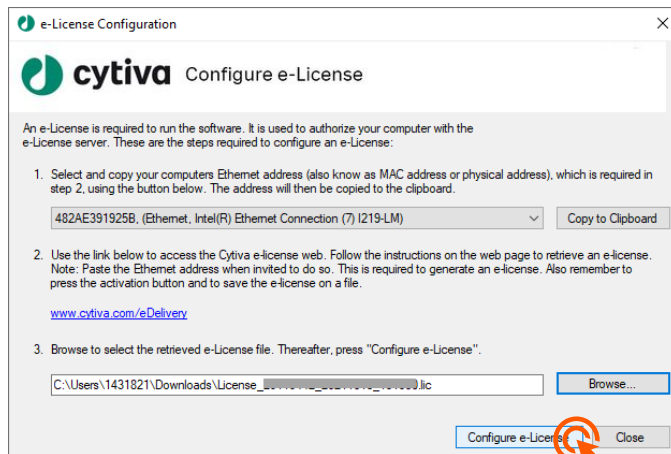


Step	Action
------	--------

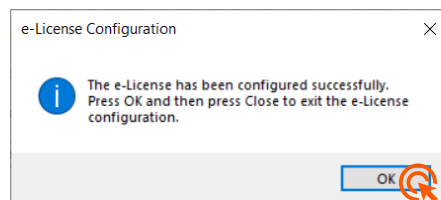
- | | |
|----|-----------------------------------|
| 14 | Browse to the e-License location. |
|----|-----------------------------------|



- | | |
|----|---------------------------------------|
| 15 | Click Open . |
| 16 | Complete the e-License configuration. |

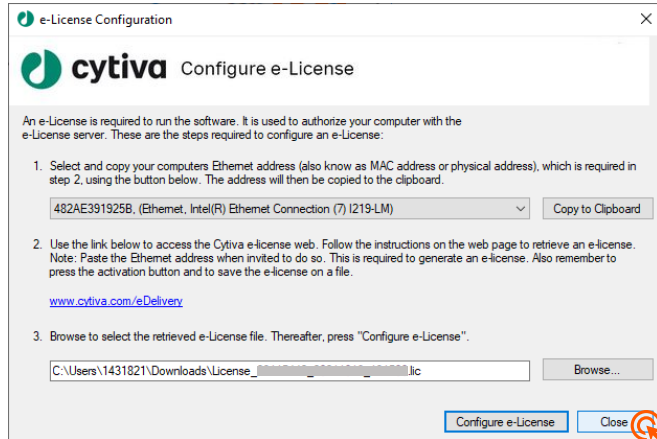


- | | |
|----|--|
| 17 | When the following dialog box appears, the configuration is successful . Click OK to continue. |
|----|--|



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

18	Close the e-License Configuration window.
----	--



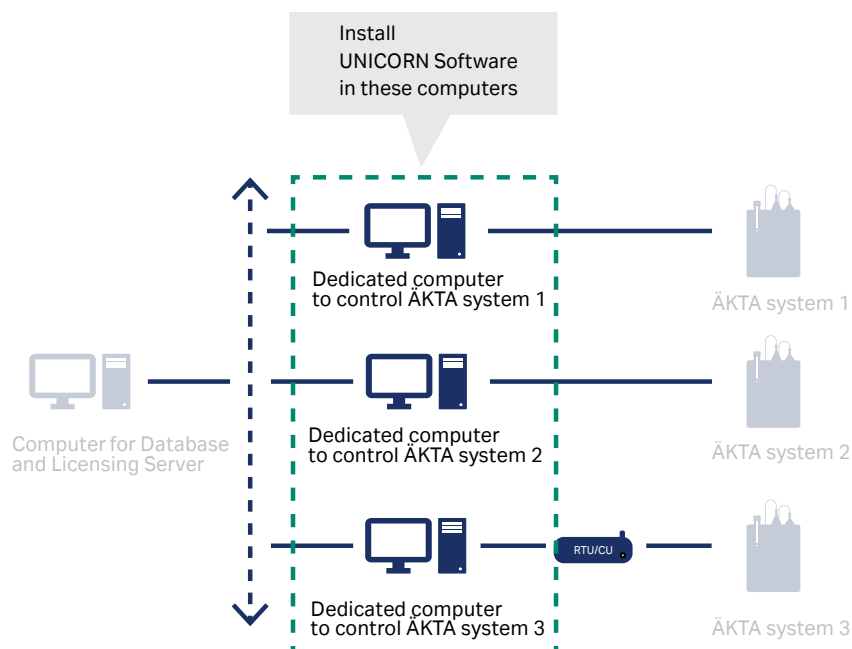
19	Go to Section 3.4 Step 4: Install the UNICORN software, on page 55 .
----	--

3.4 Step 4: Install the UNICORN software



IMPORTANT

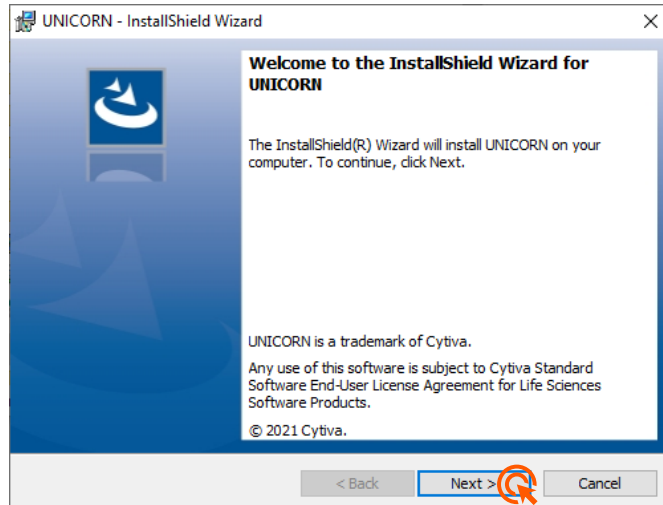
Perform the following steps in all the dedicated computers controlling your ÄKTA systems, see the illustration below.



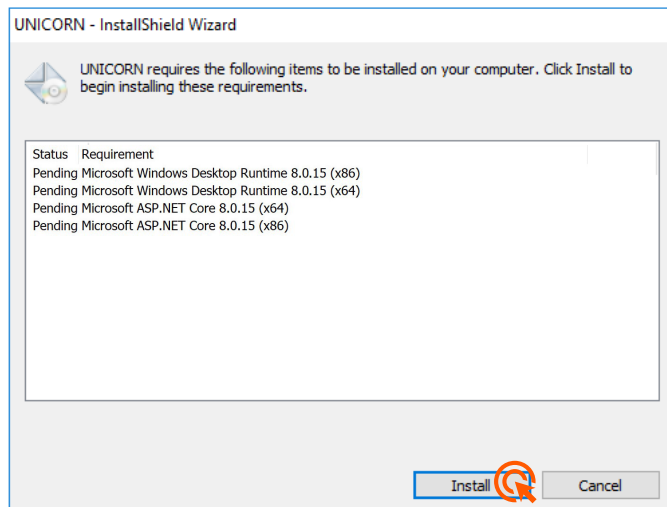
Step	Action
1	Locate and double-click the UNICORN_7.x.x.xxx_eDelivery_Portal folder.
2	Double-click Setup.exe . If the Open File - Security Warning window appears, click Run .

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

3	Click Next .
---	---------------------

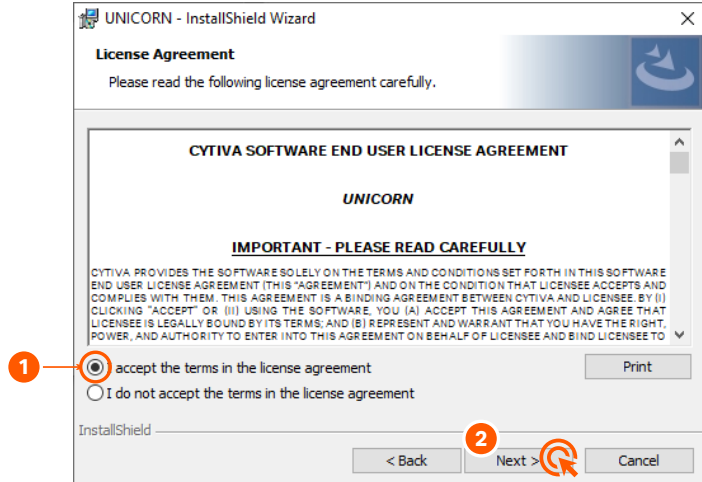


4 If your computer is **missing any software requirements**, the installer will install them. The computer can **restart** several times. After the restart, double-click the **Setup.exe** again, see **step 5**.

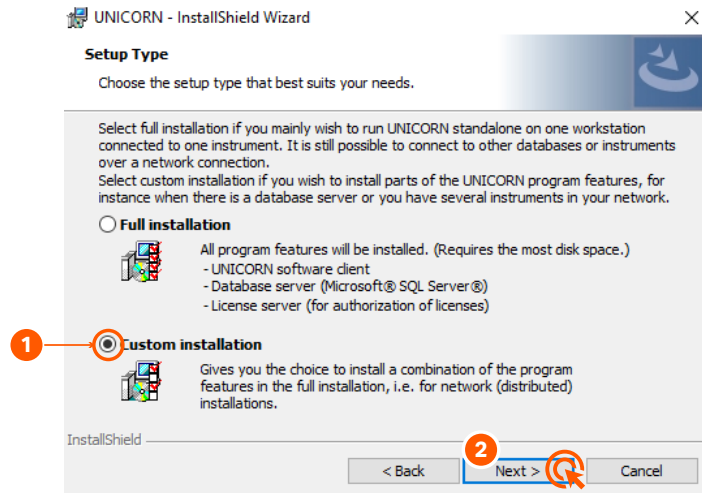


Step	Action
------	--------

- | | |
|---|--|
| 5 | Accept the license agreement and click Next . |
|---|--|

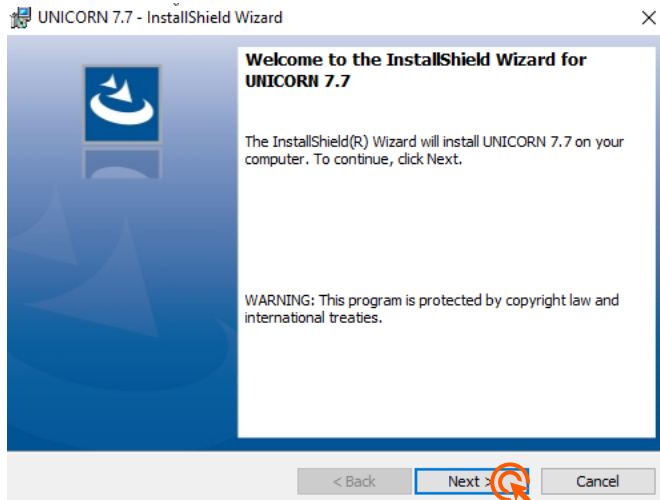


- | | |
|---|--|
| 6 | Select the installation type and click Next . |
|---|--|

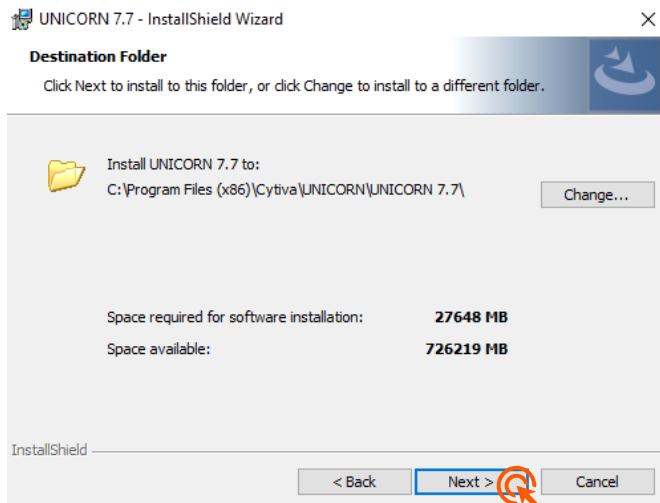


Step	Action
------	--------

7	Click Next .
---	---------------------

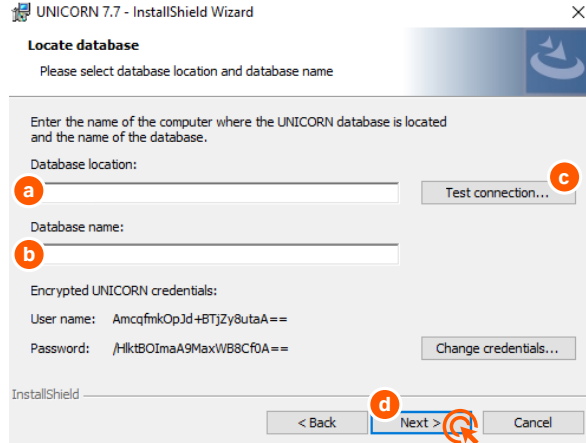


8	Click Next .
---	---------------------



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

9	Follow the steps in the illustration below.
---	---



- Enter the name of the computer where the database is installed.
- Enter the name of the database.
- Click **Test connection** to check if the database is reachable.
Result: A pop-up window appears to confirm if the connection is successful. Click **OK** on the pop-up window.
- Click **Next** to continue.

10

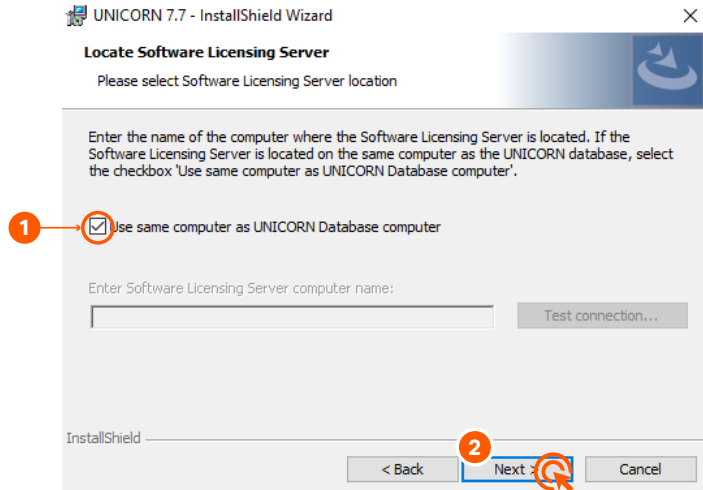


IMPORTANT

If you have chosen to install the **Software Licensing Server** on a different computer, enter the licensing server name instead of selecting **Use same computer as UNICORN Database computer**.

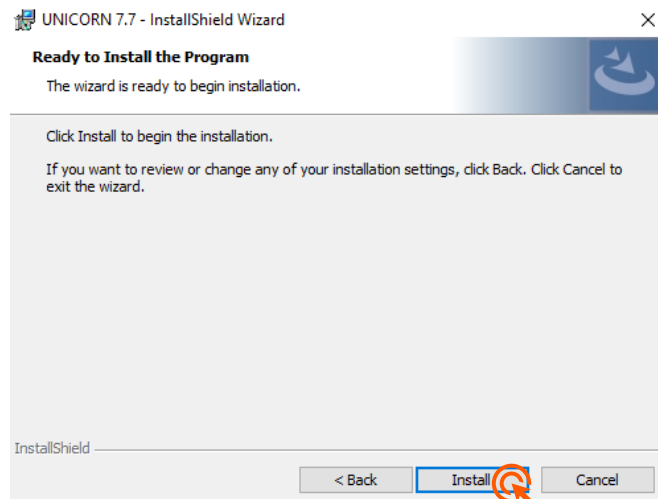
Step	Action
------	--------

Follow the steps in the illustration below.



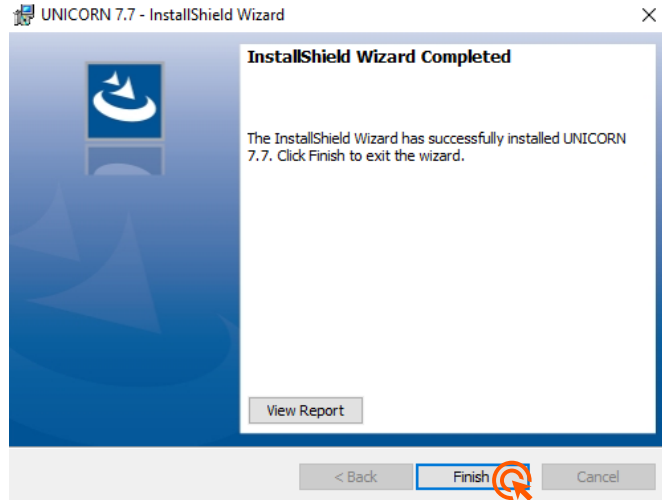
11

Click **Install**.

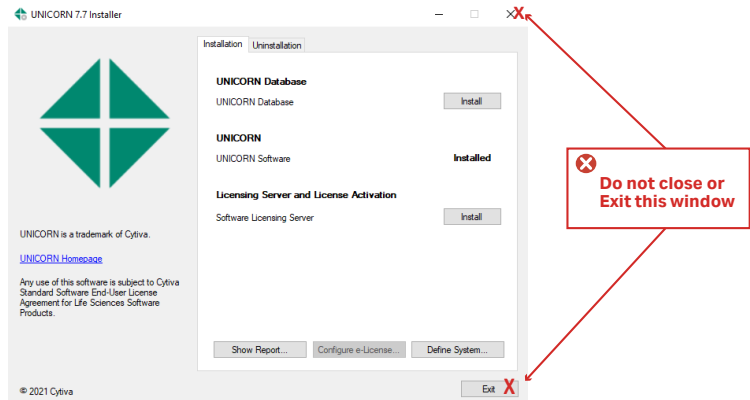


Step	Action
------	--------

12	Click Finish .
----	-----------------------



The following window confirms that the **UNICORN Software** is installed successfully.

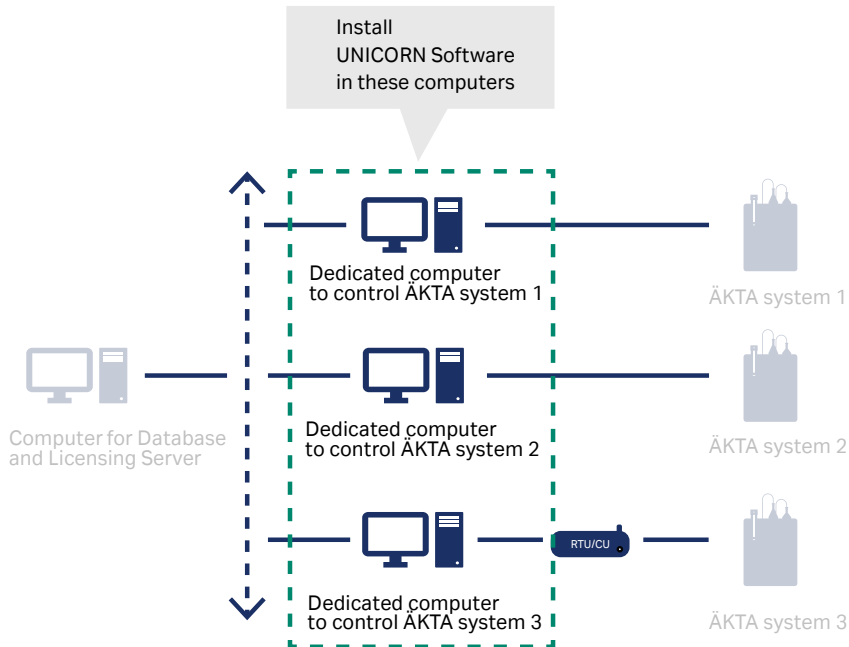


13	Go to Section 3.5 Step 5: Configure the IP, on page 62 .
----	--

3.5 Step 5: Configure the IP



IMPORTANT
Perform the following steps in all the dedicated computers controlling your ÄKTA systems, see the illustration below.

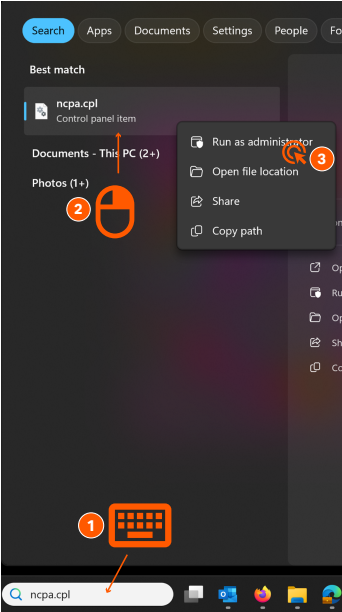


Step	Action
1	Make sure that your dedicated computer is connected to the Internet and to the ÄKTA system.
2	Click the Windows search icon.



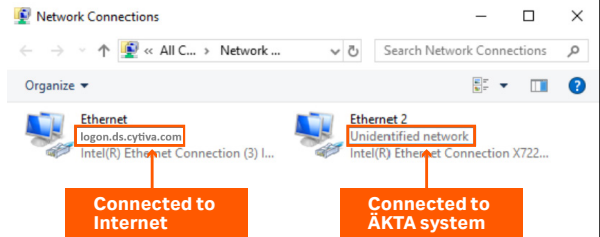
Step	Action
------	--------

3	Type <i>ncpa.cpl</i> in the search box , and then press Enter .
---	--



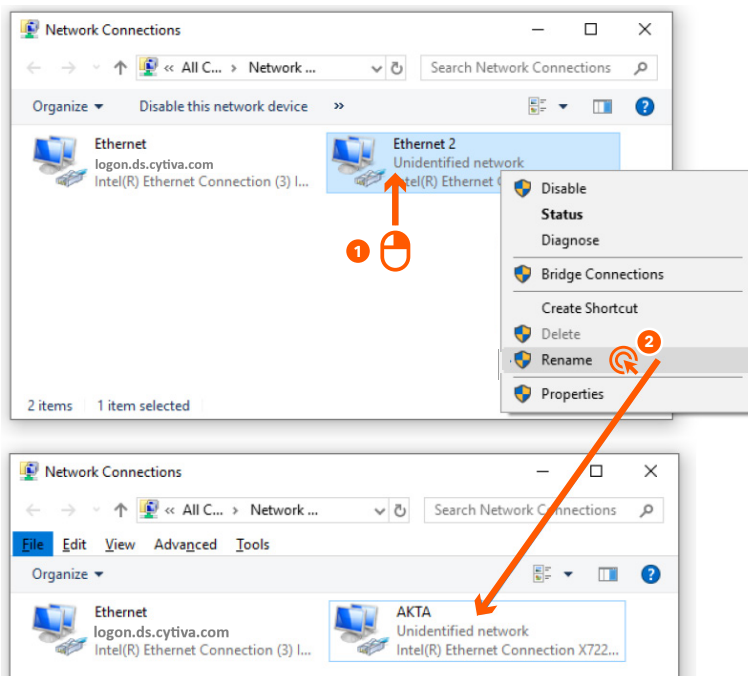
IMPORTANT

If you have two Ethernet cards, one Ethernet card is connected to the Internet (**Ethernet in the illustration below**) and the other (**Ethernet 2 in the illustration below**) is connected to the ÄKTA system. Generally, the Ethernet card connected to the Internet has the company domain name. For example, in the illustration below, it is **logon.ds.ge.com**. The Ethernet card connected to the ÄKTA system generally states **Unidentified network** as seen in the illustration below.

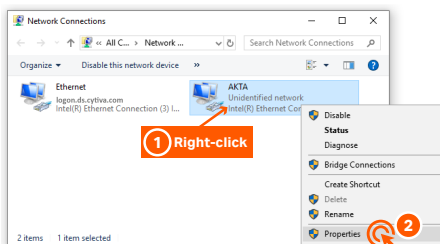


Step	Action
------	--------

- | | |
|---|--|
| 4 | Rename the Ethernet 2 connection to AKTA . |
|---|--|

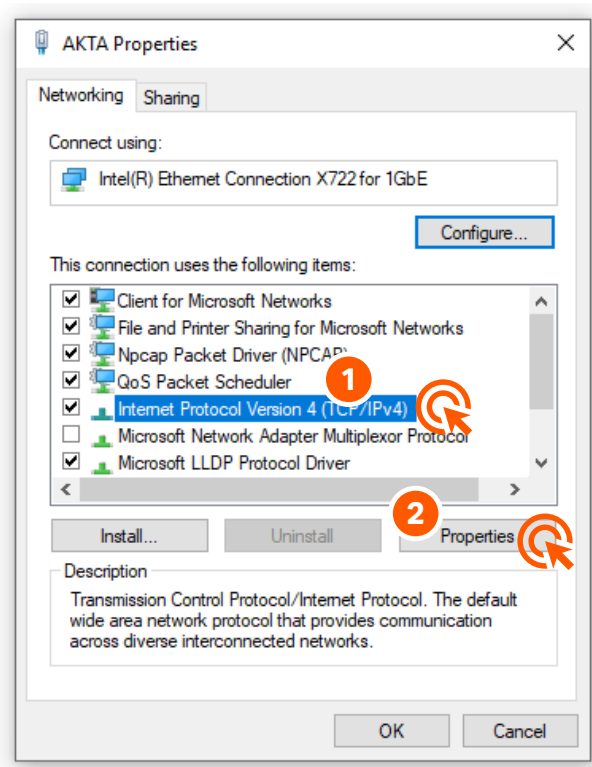


- | | |
|---|---|
| 5 | Right-click the AKTA network connection and select Properties . |
|---|---|



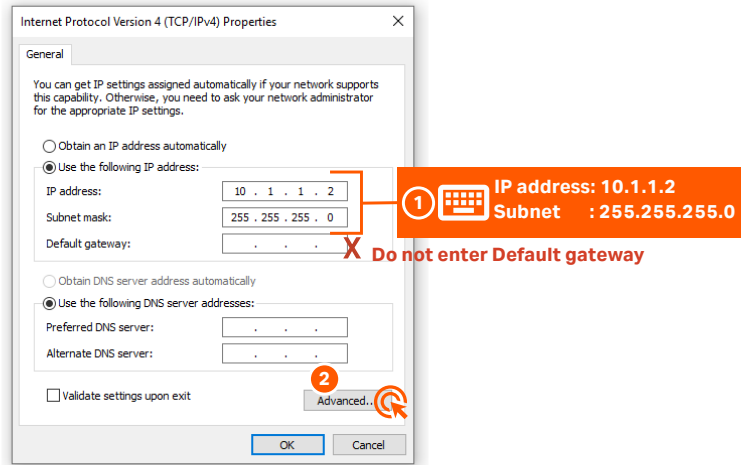
Step	Action
------	--------

6	Select Internet Protocol Version 4 .
---	---

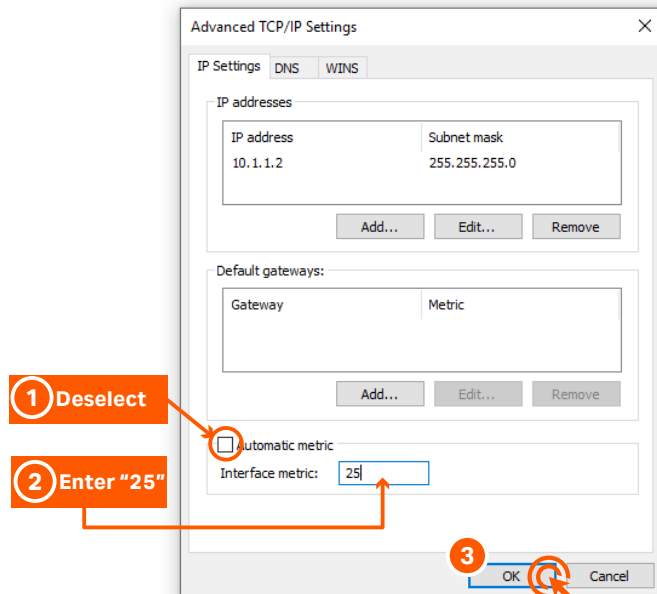


Step	Action
------	--------

- | | |
|---|--|
| 7 | Enter the IP address and the Subnet mask as shown in the illustration below. |
|---|--|



- | | |
|---|--|
| 8 | Apply the advanced TCP/IP settings as shown in the illustration below. |
|---|--|



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

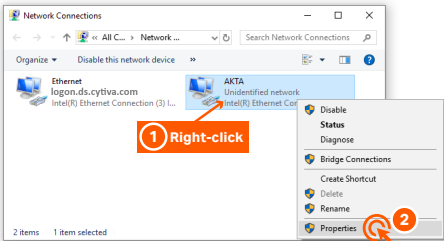
9	Click OK or Close in all the open dialogs.
---	--



IMPORTANT

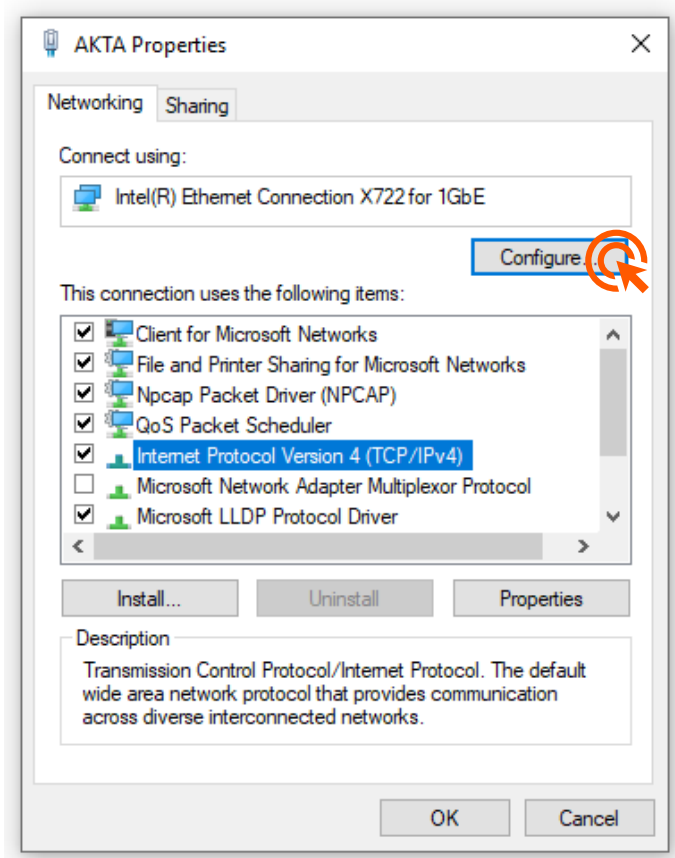
- If the system to be connected is a CU-based system, for example, ÄKTApromess™, ÄKTA ready™, ÄKTA readyflux™, etc., perform steps 10 to 12.
- If the system to be connected is not a CU-based system, for example, ÄKTA avant, ÄKTA pure™, ÄKTA go™, ÄKTA pilot™ 600, ÄKTA process™, etc., skip steps 10 to 12. The **Speed & Duplex** property for these systems is set to **Auto Negotiation** by default.

10 Access the **Properties** window of the ÄKTA network connection.



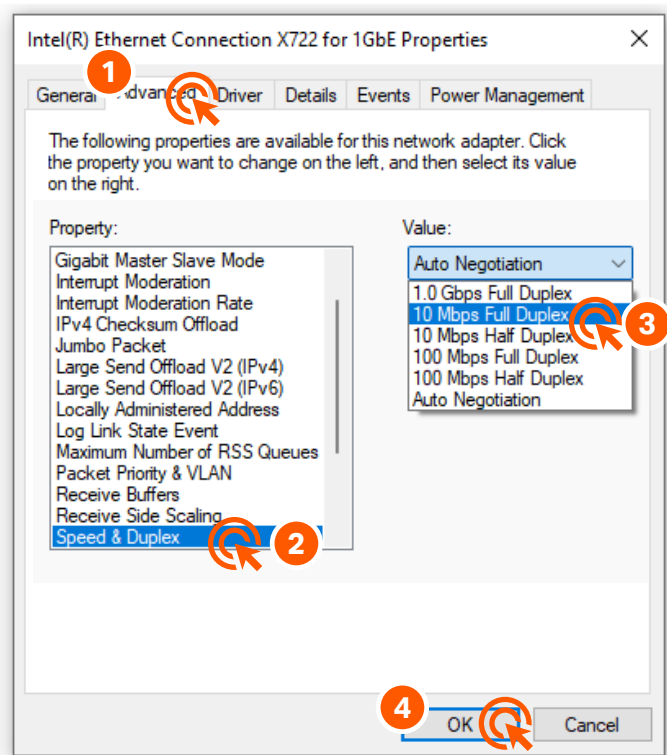
Step	Action
------	--------

11	Click Configure...
----	---------------------------



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

- | | |
|----|--|
| 12 | Set the Speed & Duplex property value to 10 Mbps Full Duplex as shown in the illustration below. |
|----|--|



- | | |
|----|---|
| 13 | Go to Section 3.6 Step 6: Define a system, on page 70 . |
|----|---|

3.6 Step 6: Define a system

See [Section 2.5 Step 5: Define a system, on page 31](#).

4 Troubleshooting

For information on troubleshooting, see the *UNICORN Administration and Technical Manual*. This document is available on the Cytiva web.

**Give feedback on this document**

Visit cytiva.com/techdocfeedback or
scan the QR code.



cytiva.com/unicorn

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

ÄKTA, ÄKTA go, ÄKTA pilot, ÄKTA pure, ÄKTAprocess, ÄKTA process, ÄKTA ready, ÄKTA readyflux, and UNICORN are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

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

© 2020-2025 Cytiva

UNICORN © 2020-2025 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

29414475 AD V:5 06/2025