



UNICORN™ Service Tool

User Manual

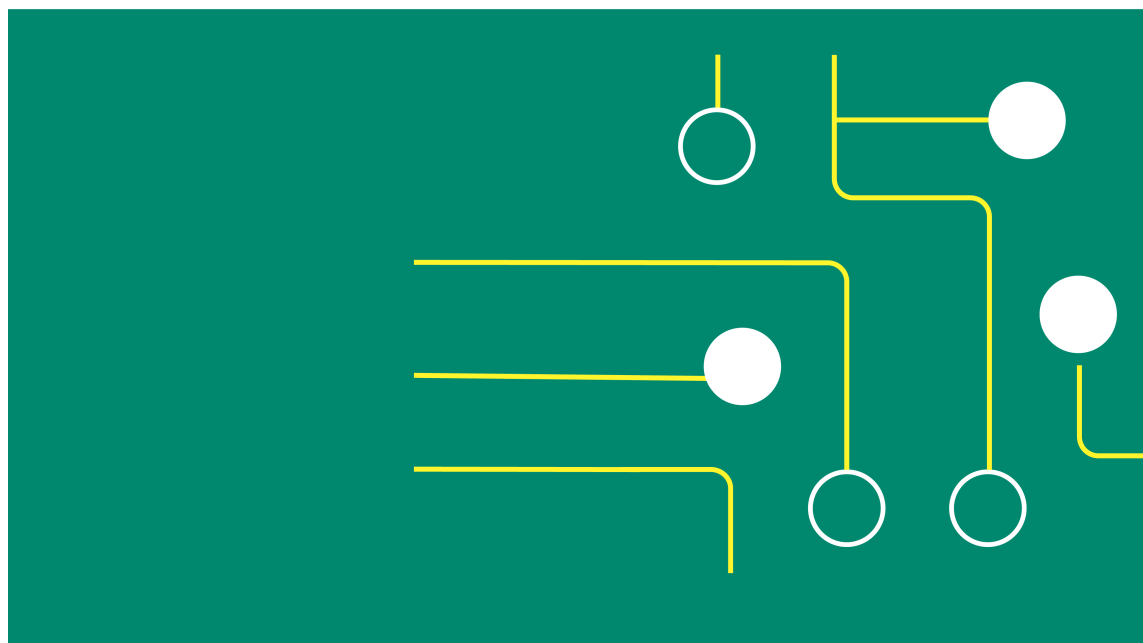


Table of Contents

1	Introduction	3
2	UNICORN Installation tab	5
3	UNICORN Connection tab	6
4	Database Server Connection tab	11
5	Optional Databases	14
6	License Server Connection tab	17
7	Active Directory Connection tab	18
8	UNICORN online tab	19
9	Create System Report tab	21
10	Process Status tab	22
11	Port check tab	24
12	Computer HW/SW check tab	29
13	Deployment tab	31
14	History Log tab	33
15	My Instruments tab	34
16	About tab	35

1 Introduction

Introduction to the UNICORN Service Tool

The UNICORN™ system control software Service Tool is used to troubleshoot and set parameters in an existing UNICORN installation. It can also be used to make sure that the operating system environment is properly set up prior to installing UNICORN.

With the UNICORN Service Tool, it is possible to do the following:

- See a list of installed UNICORN components
- Configure IP addresses of the *Real-Time Unit* (RTU)
- Test connection to a database server
- Test connection to a license server
- Manage processes needed by UNICORN
- Test communication ports
- Check if the computer and operating system settings fulfills the UNICORN specifications
- See a log of all events registered by the UNICORN Service Tool
- Update RTU operating system and software

Typographical conventions

Menu commands, field names and other text items from the software are quoted exactly as they appear on the screen, in a bold italic typeface:

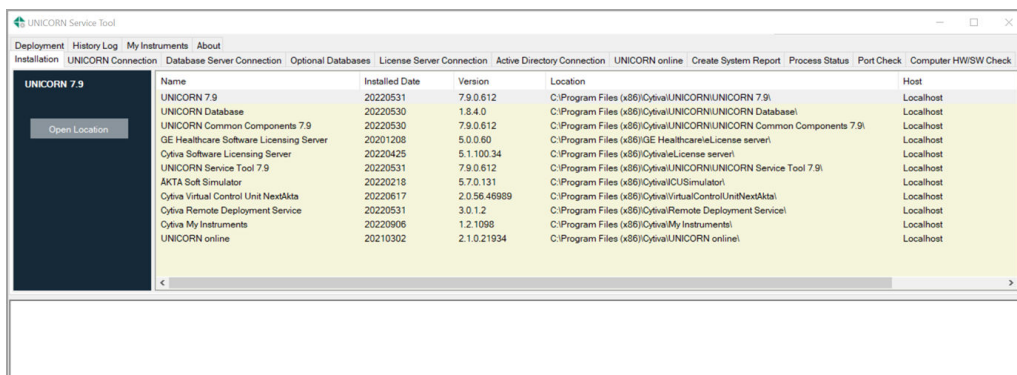
Example: ***Port check*** tab

Start UNICORN Service Tool

Step	Action
1	Click the Windows <i>Start</i> button.
2	Type UNICORN Service Tool in the <i>Start</i> menu search field.
3	Click the <i>UNICORN Service Tool</i> item that is displayed as the search result to start the program.

The UNICORN Service Tool user interface

The different functions of the UNICORN Service Tool are accessed by clicking the tabs of the UNICORN Service Tool window. There is a status pane in the lower part of the window.



The screenshot shows the UNICORN Service Tool window. The 'UNICORN 7.9' tab is selected in the left sidebar. The main area displays a table of installed components.

Name	Installed Date	Version	Location	Host
UNICORN 7.9	20220531	7.9.0.612	C:\Program Files (x86)\Cyteva\UNICORN\UNICORN 7.9\	Localhost
UNICORN Database	20220530	1.8.4.0	C:\Program Files (x86)\Cyteva\UNICORN\UNICORN Database\	Localhost
UNICORN Common Components 7.9	20220530	7.9.0.612	C:\Program Files (x86)\Cyteva\UNICORN\UNICORN Common Components 7.9\	Localhost
GE Healthcare Software Licensing Server	20201208	5.0.0.60	C:\Program Files (x86)\GE Healthcare\License server\	Localhost
Cytiva Software Licensing Server	20220425	5.1.100.34	C:\Program Files (x86)\Cytiva\License server\	Localhost
UNICORN Service Tool 7.9	20220531	7.9.0.612	C:\Program Files (x86)\Cytiva\UNICORN\UNICORN Service Tool 7.9\	Localhost
AKTA Soft Simulator	20220218	5.7.0.131	C:\Program Files (x86)\Cytiva\ICUSimulator\	Localhost
Cytiva Virtual Control Unit NextAkta	20220617	2.0.56.46989	C:\Program Files (x86)\Cytiva\VirtualControlUnitNextAkta\	Localhost
Cytiva Remote Deployment Service	20220531	3.0.1.2	C:\Program Files (x86)\Cytiva\Remote Deployment Service\	Localhost
Cytiva My Instruments	20220906	1.2.1098	C:\Program Files (x86)\Cytiva\My Instruments\	Localhost
UNICORN online	20210302	2.1.0.21934	C:\Program Files (x86)\Cytiva\UNICORN online\	Localhost

Tab
Installation
UNICORN Connection
Database Server Connection
Optional Databases
License Server Connection
Active Directory Connection
UNICORN online
Create System Report
Process Status
Port check
Computer HW/SW check
Deployment
History Log
My Instruments¹
About

¹ Only available if My Instruments 1.2 is installed.

2 UNICORN Installation tab

Introduction

Click the **UNICORN Installation** tab to see a list of installed UNICORN components. If the Service Tool is connected to an RTU, some installations can be followed by their Host IP addresses.

Copy the list of installed UNICORN components

Step	Action
1	Right-click an item in the list.
2	Click: • Copy to copy the selected rows to the Clipboard. - or • Copy Table to copy the entire list to the Clipboard.

Open an installation folder of a UNICORN component

The following steps are applicable for local installations only.

Step	Action
1	Select an item in the list.
2	Click Open Location . <i>Result:</i> The installation folder of the selected item opens in Windows Explorer.

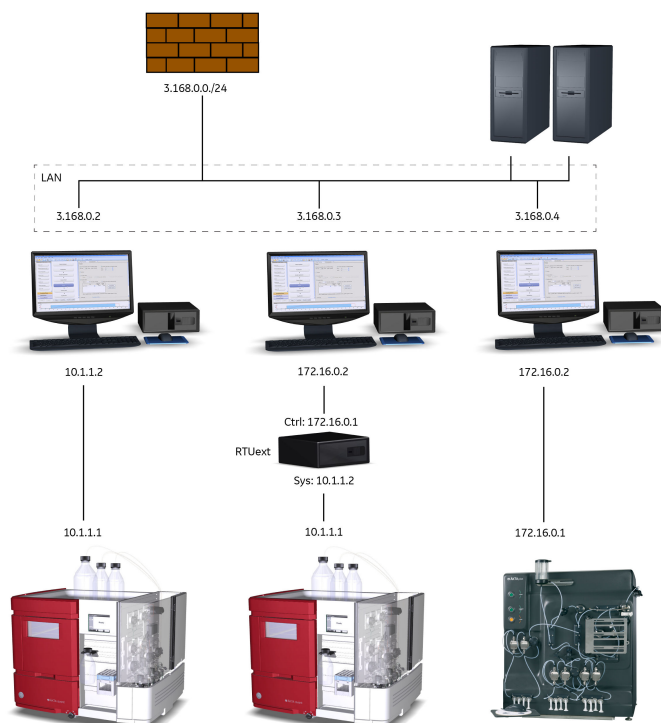
3 UNICORN Connection tab

Introduction

Select the **UNICORN Connection** tab to do the following:

- Configure IP addresses of RTU
- Search for connected instruments
- Update RTU operating system and software

The following illustration is an example network and the basic principles can be followed when configuring IP addresses:



RTU settings

The RTU has two network adapters. One connects the UNICORN control PC and the other connects the system. The Service Tool can be used to configure the IP addresses that the RTU can have on these adapters.

The **Network adapter(s)** list under **Real-Time Unit Settings** contains the connected network adapters to detect the RTUs. When a detected RTU is selected from this list, the following appear:

- A set of valid IP addresses to be used for UNICORN control PC communication appears in the **Real-Time Unit IP address - Ctrl** list. The selected IP address must be on the same subnet as the network adapter on the UNICORN control PC used for system communication.
- A set of valid IP addresses to be used for system communication appears in the **Real-Time Unit IP address - Sys** list. The selected IP address must be on the same subnet as the IP address of the system. The list appears disabled if only one IP address is available. The set of valid IP addresses are listed in the Service Tool configuration file.

- Note:**
- *The IP address of the RTU must be changed before changing the IP address of the local network adapter.*
 - *If the IP address for UNICORN control PC is changed, the connection to the RTU is lost until the local network adapter is updated to a matching IP address.*

Step	Action
1	Select an IP address for the UNICORN control PC communication from the RTU IP address - Ctrl list.
2	Optionally, select an IP address for system communication from the RTU IP address - Sys list.
3	Click Apply to configure the RTU.
	Tip: • <i>If a system is connected prior to the IP address change, use the Edit dialog to update the field RTU IP address for the system.</i>

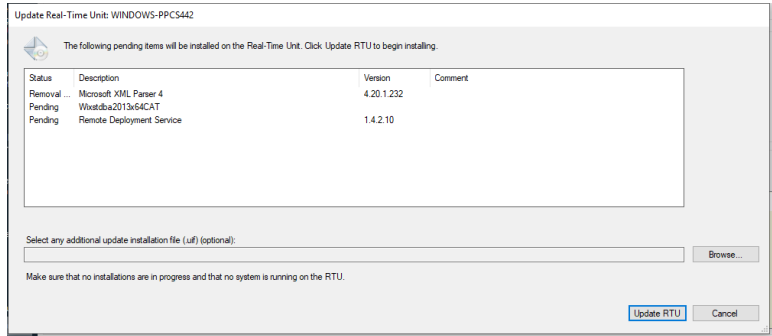
Update RTU operating system and software

Do the following steps to update to new versions of the RTU operating system, software, and service packs:

Step	Action
1	Select the Connection tab in UNICORN Service Tool.

Step Action

- Click **Update** to see the software available for update. All files will be updated in the order as shown in the list. If there are no available updates to install, go directly to the next step.

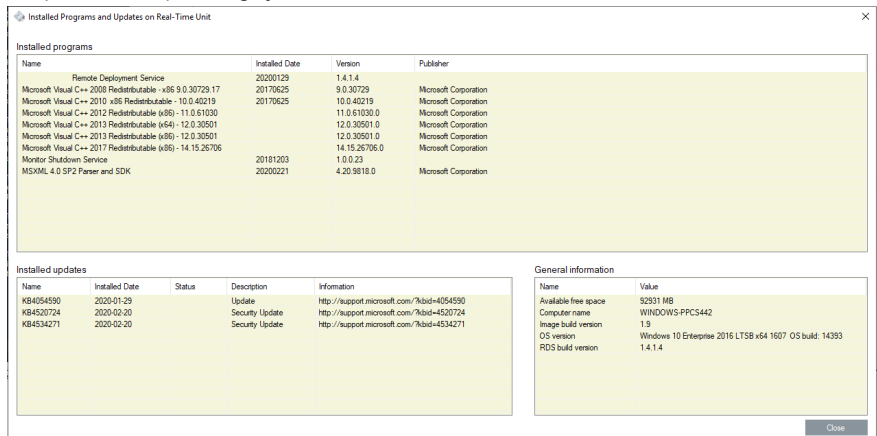


- Browse to the location of the additional downloaded software and service packs to install.
- Click **Update RTU** to start the update of the listed software and service packs. You can see the progress of the update in the log window.

View details

Click **Details** to see the software that is already installed.

After that you can click **Update** to start installation of available updates as described in *Update RTU operating system and software*.



Reset update

If something does not turn out as expected during software update you can reset some of the installed software to the point as before the update. Reset can only be applied to UNICORN control software, for example, LS Gateway, Virtual Control Unit, and Virtual Control Unit NextAkta.

Click **Reset** to make a rollback to reset the software update.

After that you can click **Update** to start installation of available updates again as described in *Update RTU operating system and software*.

Search for a connected instrument

It is possible to search for connected instruments using the Service Tool. Instruments connected directly to the UNICORN control PC, via a locally connected RTU, or connected on the same subnet on a network can be found. The search is performed by clicking the **Search** button. All instruments found are listed with the following information:

Instrument Serial Number	The serial number of the instrument
Status	Connected/Not connected to an RTU
IP Address	IP address of the instrument
Found on network	The IP address of the local network adapter that the instrument is found on
Used by UNICORN from	The IP address of the connected computer

Step	Action
1	Click Search to search for instruments connected to the selected network adapter. <i>Result:</i> Connected instruments are listed in the Found instruments list.

Copy the list of connected instruments

Step	Action
1	Right-click an item in the list.
2	Click: a. Copy to copy the selected rows to the Clipboard. b. Copy Table to copy the entire list to the Clipboard.

Configure instrument connection timeout

If the connected instrument keeps losing connection, configure the timeout period.

Note: Use higher timeout values for troubleshooting only. The default value is the only recommended setting. Higher timeout values can result in delayed display of events and response to commands.

Note: Change this setting only if all other issues, such as incorrectly setup firewall, have been ruled out.

Note: The settings are only valid on the UNICORN control PC.

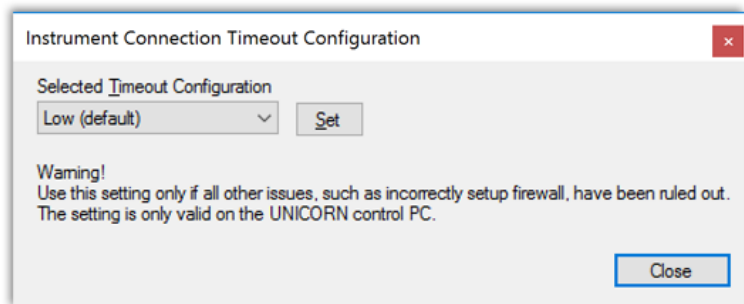
Value	Timeout period (s)
Low (default)	2
Medium	15
High	30

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

1	In the left pane of the Unicorn Connection tab, click Configure .
---	---

Result:

The **Instrument Connection Timeout Configuration** dialog box opens.



2	In the Selected Timeout Configuration list, select Medium , then click Set .
---	---

3	Restart the UNICORN control PC.
---	---------------------------------

4	If the Medium setting does not resolve the issue, select the High setting.
---	--

Tip: If a higher timeout value resolves the issue it can be an indicator for an unstable network.

4 Database Server Connection tab

Introduction

Click the **Database Server Connection** tab to search for UNICORN databases and configure UNICORN.

Search for UNICORN databases

Click **Search for UNICORN databases** to search for UNICORN databases on the network.

Result: Found databases are listed in the box below the button. The lists **Select database location** and **Select database name** are populated with found database locations and database names.

Note: *In order to find databases from UNICORN versions 7.6, 7.7, 7.8, 7.9, and 7.10, the latest UNICORN Service Packs must be installed on the computers holding those databases.*

Databases from UNICORN 7.5 and earlier versions cannot be found.

Test database connection

Click **Test Database Connection** to test the connection to the selected database.

Result: If the database connection is successful, the icon to the right of the **Test Database Connection** button becomes green and the lower status pane is updated with the result of the test.

Configure UNICORN

Step	Action
1	Select a database location in the Select database location list or enter the computer name where the database is located.
2	Select a database name in the Select database name list or enter a database name.
3	a. Turn on or turn off encrypted connection to the database by selecting or clearing the Encrypt connection check box.

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

- | | |
|--|--|
| | <p>b. If the database server does not have a trusted certificate issued by a Certificate Authority, select the Trust server certificate check box. This check box is selected by default.</p> |
|--|--|

Note:

The **Trust server certificate** check box is only available if the **Encrypt connection** check box is selected.

- | | |
|---|--|
| 4 | Click Configure UNICORN to configure UNICORN to use the selected database with the selected security options. |
|---|--|

Configure database user name and password

Configure the database login credentials used by UNICORN. The credentials can be entered in plain text which are automatically encrypted.

Note: *Encrypted user name and password are typically provided by the database administrator.*

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

- | | |
|---|--|
| 1 | Click Edit to open the Configure UNICORN Database credentials dialog box. |
| 2 | <ul style="list-style-type: none"> Enter a new user name in the User name box. Enter a new password in the Password box. |

Note:

The **Encrypted credentials** will be populated automatically.

Or

Enter the encrypted credentials in the relative boxes.

- | | |
|---|---|
| 3 | Click Apply to configure UNICORN to use the new encrypted database password. |
|---|---|

Note: *Do not use the **Configure UNICORN Database credentials** dialog box unless it is certain that the database password has been changed by the database administrator. Change only the password that UNICORN uses to access the database, not the actual database password.*

5 Optional Databases

Introduction

Optional Databases tab is used to enable the function that allows the user to select which database to log on to when starting UNICORN. The list of databases that can be selected when starting UNICORN is defined in this tab and UNICORN is configured accordingly.

Enable Optional Databases

Step	Action
1	Enable or disable Optional Databases by selecting or deselecting the Enable Optional Databases check box.
2	Click the Configure UNICORN button to validate and apply the change.

Test connection

Click **Test Connection** to test the connection to the selected database server.

Result:

If the database connection is successful, the little white circle next to the **Test Connection** button becomes green and the status pane next to it is updated with the result of the test.

Define UNICORN databases

The list of databases that can be selected when starting UNICORN can be defined in the following ways:

- Automatic search for UNICORN Databases
- Default manual definition of UNICORN Databases
- Advanced manual definition of UNICORN Databases

Automatic search for UNICORN databases

Step	Action
1	Click Search for UNICORN Databases to search for UNICORN Databases on the network.
2	In the Select Databases To Add window, select databases and click Add .

Note: In order to find databases from UNICORN versions 7.6, 7.7, 7.8, 7.9, and 7.10, the latest UNICORN Service Packs must be installed on the computers holding those databases.

Databases from UNICORN 7.5 and earlier versions cannot be found.

Default manual definition of UNICORN databases

Step	Action
1	Enter an alias in the Alias text box (optional).
2	Enter a server location in the Server text box, for example: <code>LOCALHOST \UNICORN</code> .
3	Click the Add button to add the database to the Optional Database List . If needed, select databases from this list and click Remove to remove them from the list.
4	Click the Configure UNICORN button to configure the UNICORN client with the added databases.

Tip:

Click **Refresh** to refresh the list.

Advanced manual definition of UNICORN databases

Step	Action
1	Select the Advanced check box.
2	Enter an alias in the Alias text box (optional).
3	Enter a server location in the Server text box, for example: <code>LOCALHOST \UNICORN</code> .
4	Enter a database name in the Database text box.
5	Enter a database user name in the User Name text box.
6	Enter the database user password in the Password text box.

Note:

The **User Name** and/or **Password** can be entered as either encrypted or in plain text.

- | | |
|---|--|
| 7 | <ul style="list-style-type: none"> Turn on or turn off encrypted connection to the database by selecting or clearing the Encrypt connection check box. |
|---|--|

Step	Action
	• If the database server does not have a trusted certificate issued by a Certificate Authority, select the Trust server certificate check box. This check box is by default selected. Note: <i>The Trust server certificate check box appears available when the Encrypt connection check box is selected.</i>
8	Click the Add button to add the database to the Optional Database List . If needed, select databases from this list and click Remove to remove them from the list.
9	Click the Configure UNICORN button to configure the UNICORN client with the added databases.
	Tip: <i>Click Refresh to refresh the list.</i>

6 License Server Connection tab

Introduction

Click **License Server Connection** to search for a license server and configure UNICORN.

Search for license server

Click **Search for Licenses** to search for licenses on the selected computer and on the localhost.

Result: Found licenses are listed in the **Found licenses** box.

Open license file location

There are two folders where the license files are stored.

- Click the upper **Open Location** button to open the folder, where the node locked license files are stored, in Windows Explorer.
- Click the lower **Open Location** button to open the folder, where the concurrent license files are stored, in Windows Explorer.

Configure UNICORN

Step	Action
1	Select a computer name in the Select or enter computer name list or enter the computer name where the license server is running.
2	Click Configure UNICORN to configure UNICORN to use the license server running on the selected computer.
Note: <i>Any node locked license on the localhost is always available in UNICORN.</i>	

7 Active Directory Connection tab

Introduction

Click the **Active Directory Connection** tab to configure the connection between UNICORN and Active Directory.

Test Active Directory connection

Step	Action
1	Select an Active Directory in the list or enter an Active Directory name.
2	Click Test Connection to test the connection to the selected Active Directory with the selected security options. <i>Result:</i> If the Active Directory connection is successful, the icon to the right of the Test Connection button becomes green.

Configure UNICORN

Step	Action
1	Turn on or turn off encrypted connection to the Active Directory by selecting or clearing the Encrypt connection check box.
2	Enter a connection port in the Port box to be used for encrypted Active Directory connection. The Default port is 636. Note: <i>The Port box is only available if the Encrypt connection check box is selected.</i>
3	Click Configure UNICORN to configure UNICORN to use the selected security options for the Active Directory connection.

8 UNICORN online tab

Introduction

UNICORN online is a web-based application that interacts with UNICORN through certain entry points. For this interaction to work, certain ports and certificates has to be configured in UNICORN. All the configurations can be done on the **UNICORN online** tab.

Enable UNICORN online application

1. Enable or disable **UNICORN online** application by selecting or deselecting the **Enable** check box.
2. Click **Configure UNICORN online** button to validate and apply the change.

Generate and configure UNICORN SSL certificates

UNICORN SSL certificate is used to encrypt the data sent between UNICORN and **UNICORN online**.

Select either of the following options:

- **Standard (Generate SSL certificates)**
- **Custom (Generate SSL certificate from issuer certificate file)**

Standard (Generate SSL certificates)

1. Select **Standard (Generate SSL certificates)**.
2. Click **Generate SSL certificate**.

Result: New certificate is generated, and UNICORN configuration is updated.

Custom (Generate SSL certificate from issuer certificate file)



IMPORTANT

To be able to generate a new certificate, you need to have an issuer certificate available.

Do the following to generate a new certificate:

Step	Action
1	Select Custom (Generate SSL certificate from issuer certificate file) .
2	Click Generate SSL certificate .
3	Browse and select the issuer certificate (PFX file format).
4	Enter password for the issuer certificate and click Create . <i>Result:</i> A new SSL/TLS certificate is created, and the new certificate information is shown under Created UNICORN SSL certificate .
5	Click Apply to use the created certificate. <i>Result:</i> The UNICORN configuration is updated.
6	Restart the computer to activate the new certificate.

Change port number

The **Ports** tab has information about all the ports used by UNICORN online. Here it possible to select which port to use if it conflicts with another application process with UNICORN on the local network (for example, used for web hosting, result viewer, authentication, and system control).

Do the following to change port number:

1. To change to the default value, go to **Port number** and doubleclick the line with the value shown in parenthesis (xxx). Or, type a desired port value.
2. Click the **Configure UNICORN online** button to validate and apply the change. The validation checks that the selected port values are not used by another application with UNICORN. It also checks that the value is within a valid range (1 – 65535).

9 Create System Report tab

Introduction

Click the **Create System Report** tab to create a system report including the current UNICORN configuration settings. This can be used to troubleshoot the UNICORN installation.

Create System Report

Step	Action
1	Type a description of the problem you are experiencing in the text box.
2	Click Create System Report. <i>Result:</i> The program starts to collect information about your current operating system and UNICORN configuration. This can take a few minutes. When the report file archive is created the Save System Report As dialog box opens. <i>Result:</i> Click Create System Report. The program starts to collect information about your current operating system and UNICORN configuration. This can take a few minutes. When the report file archive is created the Save System Report As dialog box opens.
3	Click Save to save the system report file archive.

10 Process Status tab

Introduction

Click the **Process Status** tab to see the list of all the UNICORN related processes. The processes can be started, stopped, registered, and unregistered in the Windows registry. If the Service Tool is connected to an EPC, some processes can be followed by their Host IP addresses.

Windows operating system settings for **service startup timeout** can be configured here.

Note: *Not all operations are available for all processes.*

Manage processes

Step	Action
1	Select a process in the list.
2	Click Start , Stop , Register or Unregister . The available buttons depend on the process and the status of the process. <i>Result:</i> The process Status changes and the lower pane is updated.

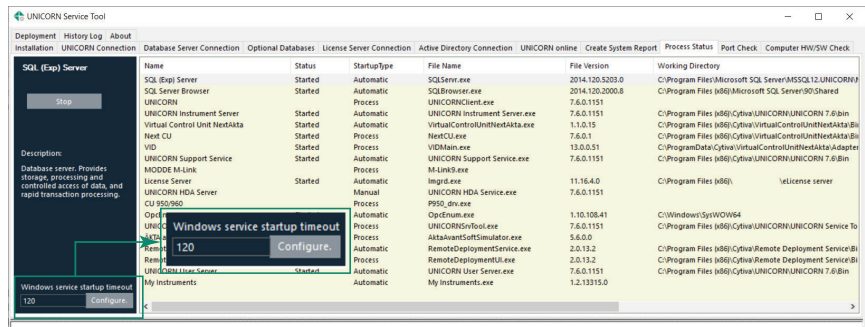
Note: *The process named **UNICORN** can not be stopped from the UNICORN Service Tool. Exit UNICORN to stop the process.*

Configure Windows services startup timeout

**IMPORTANT**
This configuration is applicable for all the services, therefore the new configuration will affect all the services.

Enter a startup timeout and click **Configure....** This is to ensure that all the services have enough time to start up before the operating system shuts them down for being too slow.

By default, the timeout is set to 30 seconds, but we recommend setting the timeout to **120 seconds**.



11 Port check tab

Introduction

The port check is intended for checking if it is possible to communicate on ports used in a UNICORN network installation. The port check verifies that computers in the network can communicate using the selected ports.

The port check can also be done locally on a single computer. That is, the port check is set up as both client and server on the local computer. This verifies that local ports can be used on the computer.

The Port check tab

Click the **Port check** tab to see a list of ports used by UNICORN and to test the connectivity of the ports.

See UNICORN Administration and Technical Manual for more information on ports used by UNICORN.

Test IP address

Before running a port check you can check that there is a connection between the client and the server computer. The port checks cannot succeed if there is no connection between the client and the server computer.

Step	Action
1	Click the Client tab.
2	Click the network adapter you want to test in the Select adapter to use list.
3	a. Enter the server computer name in the Name of computer acting server box or select a server from the list and/or b. enter the IP address in the IP address of computer acting server box.
4	Click Ping to test the IP address. <i>Result:</i> The status pane is updated with the result of the test. <i>Result:</i> Click Ping to test the IP address. The status pane is updated with the result of the test.

Step	Action
5	If the IP address test is successful, the icon to the right of the IP address of computer acting server box becomes green and the lower status pane is updated with the result of the test. If the IP address test is unsuccessful you must make sure that there is a connection between the client and the server computer.

Test client and server ports

Follow the instructions to run a port check. If the server and client are on different computers the port check must be run on both computers simultaneously.

Step	Action
1	Click the Server tab.
2	Click the network adapter connected to the client computer in the Select adapter to use list.
3	Select the ports to test by selecting the check boxes for the server ports in the left pane.
4	Click Start port check .
5	Click the Client tab.
6	Click the network adapter connected to the server computer in the Select adapter to use list.
7	Select the same ports that were selected on the Server tab by selecting the check boxes in the left pane.
8	Click Start port check. <i>Result:</i> The Status column is updated according to connection status and the lower pane is updated with the result of the test. <i>Result:</i> Click Start port check. The Status column is updated according to connection status and the lower pane is updated with the result of the test. Note: <i>If the software asks you to stop a running service when performing a port check, stop the service in the Process Status tab and then click Start port check again.</i>

Port check between computers

Different computers can have different roles in the network. For example the UNICORN Software can be installed on one computer, the Database server on another and License Server on a third computer.

In the following example the Database Server is installed on one computer and the UNICORN Software is installed on a different computer. In order to successfully complete a port check you must first start a port check on the server computer and then start a port check on the client computer.

Stage	Description
1	Start a port check on the server computer.
2	Start a port check on the client computer.
3	The port check is successful if the port check on the client computer is successful.

Start port check on the server computer

Step	Action
1	Start the UNICORN Service Tool on the Database server computer.
2	Click the Port check tab.
3	Click the Server tab.
4	Click the network adapter connected to the client computer in the Select adapter to use list. <i>Result:</i> The Name of computer acting as server and IP address of computer acting as server boxes are populated with the Database server information. Copy this information for use in the client computer. <i>Result:</i> Click the network adapter connected to the client computer in the Select adapter to use list. The Name of computer acting as server and IP address of computer acting as server boxes are populated with the Database server information. Copy this information for use in the client computer.
5	Select the SQL Database check box in the left list. <i>Result:</i> The right-hand list is populated with information on the ports and process needed for the SQL database. <i>Result:</i> Select the SQL Database check box in the left list.

Step	Action
	The right-hand list is populated with information on the ports and process needed for the SQL database .
6	Click Start port check .
7	Wait for the processes to start and the Status change to Server → OK

Start port check on the client computer

Step	Action
1	Start the UNICORN Service Tool on the UNICORN Software computer.
2	Click the Port check tab.
3	Click the Client tab.
4	Click the network adapter connected to the server computer in the Select adapter to use list.
5	a. Enter the server computer name in the Name of computer acting server field or select a server from the list and/or b. enter the IP address in the IP address of computer acting server. Note: <i>Copy the computer name and IP address from the server computer.</i>
6	Click Ping to test the IP address. Note: <i>The port checks only succeed if there is a connection between the client computer and the server computer.</i> Result: The status pane is updated with the result of the test. Continue to the next step if the ping test result is successful.
7	Select SQL database in the left-hand list. Result: The right-hand list is populated with information on the ports used by the SQL database. Result: Select SQL database in the left-hand list. The right-hand list is populated with information on the ports used by the SQL database.

Step	Action
8	Click Start port check .
9	Wait for the processes to start and for the Status to change to Client → OK . Note: <i>If the port is not available the client side will show Client → Error. If, during the the port check, the port becomes available the Status changes to Client → OK.</i>
10	Click Stop to stop the port check.
11	Click Stop in the Port check tab on the server computer to stop the port check.

12 Computer HW/SW check tab

Introduction

Click the **Computer HW/SW check** tab to see lists of operating system and computer hardware recommendations and requirements. There are three different lists, one list for each installation type. The full UNICORN installation requirements are checked by selecting all three lists.

- UNICORN Software
- Database Server
- License Server

Check installation recommendations

Select the check box for an installation type.

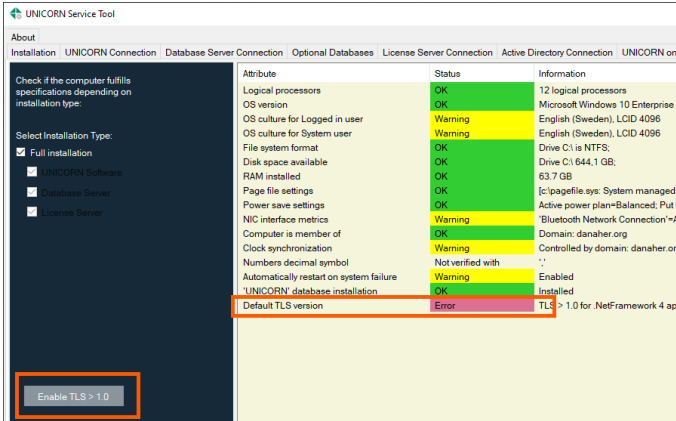
Note: Select **Full installation** to check all three installation lists.

Result: The right-hand list is populated with the recommendations for the selected installation type. The UNICORN Service Tool checks the current computer specifications against the requirements and advice user action if needed.



IMPORTANT

If you receive **Default TLS version** error, click **Enable TLS > 1.0** on the left.



Attribute	Status	Information
Logical processors	OK	12 logical processors
OS version	OK	Microsoft Windows 10 Enterprise
OS culture for Logged in user	Warning	English (Sweden), LCID 4096
OS culture for System user	Warning	English (Sweden), LCID 4096
File system format	OK	Drive C: is NTFS.
Disk space available	OK	Drive C: 644.1 GB.
RAM installed	OK	63.7 GB
Page file settings	OK	System managed
Power save settings	OK	Active power plan: Balanced
NIC interface metrics	Warning	'Bluetooth Network Connection' #2
Computer is member of	OK	Domain: danaher.org
Clock synchronization	Warning	Controlled by domain: danaher.org
Numbers decimal symbol	Not verified with	''
Automatically restart on system failure	Warning	Enabled
UNICORN database installation	OK	Installed
Default TLS version	Error	TLS > 1.0 for .NetFramework 4 ap

List of recommendations and specifications

The **Status** column for each item shows **OK** if the recommendation is fulfilled. If a specification is not fulfilled the **Status** column shows **Warning** or **Error**. Recommended user action is showed in the **Specification** column.

Note: *The list is updated when the installation type check box is selected. Clear and select the check box to update the list.*

13 Deployment tab

Introduction

The **Deployment** tab displays related information regarding the available hosts within the network using UNICORN 7.4 or later releases.

Export selected hosts

Perform these steps if you want to save information related to some hosts and later import them to another UNICORN Service Tool.

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

- | | |
|---|---|
| 1 | Select multiple hosts (Ctrl + left-click). |
| 2 | Right-click and select Export selected hosts . |
| 3 | Select a location for the file to be saved and then click Save . |

Add known host

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

- | | |
|---|--|
| 1 | Right-click and select Add known host . |
| 2 | Enter Host name / IP address and click Add .

<i>Result:</i>
The host is added to the list. |

Import known hosts

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

- | | |
|---|---|
| 1 | Right-click and select Import known hosts . |
| 2 | Locate the file created in Export selected hosts, on page 31 , and click Open .

<i>Result:</i>
The hosts listed in the file are added in the list of hosts in the Deployment tab. |

Remove selected hosts

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

1	Right-click a host and select Remove selected hosts .
---	--

Result:

The selected host is removed from the list.

Note:

Hosts that are online will reappear.

14 History Log tab

Introduction

Click the **History Log** tab to see all events registered by the UNICORN Service Tool.

Use the History Log filter

Use the History Log filter to easier find UNICORN Service Tool events.

Step	Action
1	Select the Use Filter check box. <i>Result:</i> Four more lists become available.
2	Click Level and select or clear the types of Level events to filter by.
3	Click Category and select or clear the types of Category events to filter by.
4	Click Start time to filter by events after a specific date and time.
5	Click End time to filter by events before a specific date and time.
6	Click outside the lists to apply the filter.

15 My Instruments tab

Introduction



IMPORTANT

The My Instruments tab is only available if My Instruments 1.2 is installed.

The My Instruments tab shows the validity of the certificate My Instruments is using. If the certificate is valid, the date of expiry of the certificate is also shown.

Update certificates

1. To generate new certificates and to update the My Instruments configuration, click ***Update certificates***.
2. Restart the computer to activate the new certificate.

16 About tab

The About tab

Click the **About** tab to see detailed information on the UNICORN Service Tool. Access this manual by clicking **Show Help**.

Note: *This manual is also accessible by pressing **F1** in any tab.*

**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.

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

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

© 2020-2025 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

29109881 AN V:14 04/2025