



UNICORN™ OPC Server 1.3

User Manual

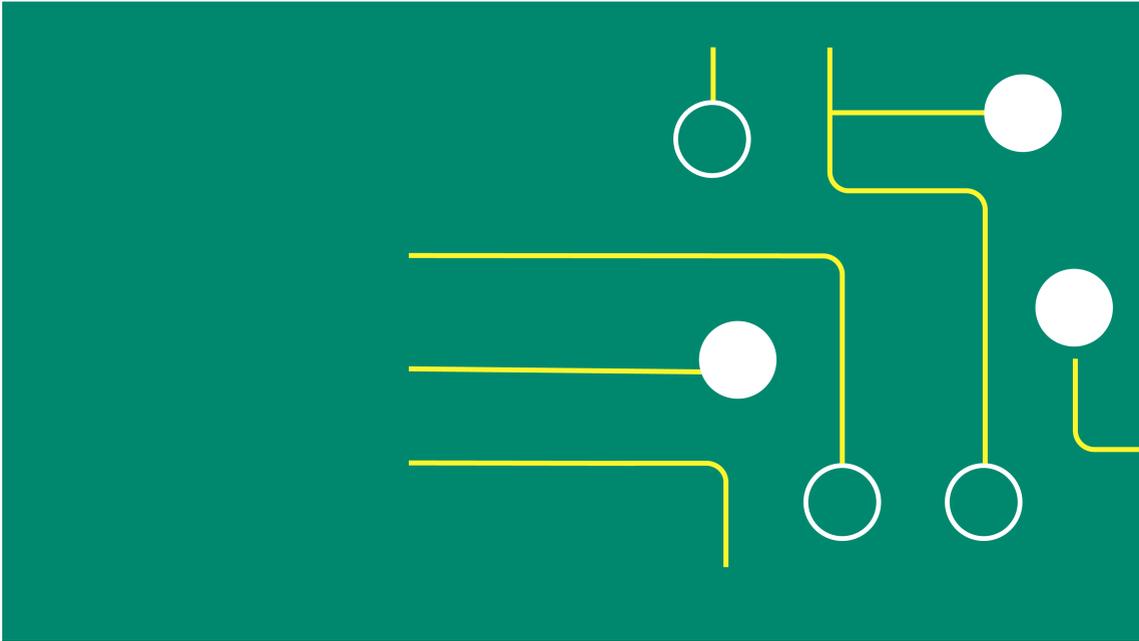


Table of Contents

1	Introduction	4
1.1	OPC Server	5
1.2	OPC UA standards and versions supported by UNICORN OPC Server	6
2	UNICORN OPC Server general settings	7
2.1	Install UNICORN OPC Server	9
2.2	UNICORN OPC Server license	11
2.3	Installation log files and server configuration files	13
2.4	Configure UNICORN OPC Server	14
2.5	Configure OPC UA Client and connect to UNICORN OPC Server	18
2.6	Trust certificate	25
2.7	Execute method/manual instruction and raise and acknowledge alarm with the OPC UA Client	28
2.7.1	<i>Connect OPC Client with Instrument</i>	29
2.7.2	<i>Execute method</i>	30
2.7.3	<i>Execute manual instruction</i>	32
2.7.4	<i>Raise and acknowledge an alarm</i>	34
2.7.5	<i>Disconnect OPC Client with Instrument</i>	38
3	Deployments	39
4	UNICORN OPC Data Access address space	42
4.1	Run data and Picture data	45
4.2	Trend data	48
4.3	Manual instructions	51
4.4	Method execution	52
4.4.1	<i>Scout Run Index</i>	53
4.4.2	<i>Run Method</i>	55
4.4.3	<i>Result Name</i>	56
4.4.4	<i>Batch ID</i>	58
4.4.5	<i>StartNotes</i>	60
4.4.6	<i>MethodNotes</i>	61
4.4.7	<i>PreCompile</i>	62
4.4.8	<i>Questions</i>	63
4.4.9	<i>Variables</i>	65
4.5	State	67
4.5.1	<i>AssignState</i>	68
4.5.2	<i>RunState</i>	70
4.5.3	<i>Command</i>	72
4.5.4	<i>InstrumentStatus</i>	74
4.5.5	<i>InstrumentStatusConnection</i>	76
4.5.6	<i>UserInControl</i>	78
4.6	Recommendations	79
5	UNICORN OPC Alarms and Events address space	80
5.1	Alarms and errors	82
5.2	Event categories	84

6	UNICORN OPC Historical Data Access address space	85
6.1	AuditTrail	87
6.2	Result file information	88
6.2.1	Curves	89
6.2.2	Documentation	90
6.2.3	Peak tables	92
6.2.4	Systems	93
6.3	UNICORN OPC Historical Data Access XML format definition	94
6.3.1	BufferPro	96
6.3.2	Calibration	97
6.3.3	EvaluationLog	98
6.3.4	Run logbook	99
6.3.5	SignatureList	100
6.3.6	SnapshotList	101
6.3.7	UsedComponents	102
6.3.8	Notes	103
6.3.9	ScoutingList	104
6.3.10	SettingsList	105
6.3.11	TextInstructions	106
6.3.12	VariableList	107
6.3.13	QuestionList	108
6.3.14	PeakTable	109
6.3.15	Unicorn raw	112
6.3.16	Columns	114
6.3.17	EvaluationProcedures	116
6.3.18	FracXY	117
6.3.19	AuditTrail	118
6.3.20	Method/Result Instrument Configuration	119
6.3.21	System	120

1 Introduction

About this chapter

This chapter contains important user information, compatibility information, and a description of the intended use of the UNICORN™ OPC Server.

In this chapter

Section		See page
1.1	OPC Server	5
1.2	OPC UA standards and versions supported by UNICORN OPC Server	6

1.1 OPC Server

The UNICORN OPC Server provides a standardized integration interface to support integration between UNICORN and other software systems, such as Laboratory Information Systems and Manufacturing Execution Systems. Customers can stream the data to a historian. OPC enables open connectivity via open standards that are created in collaboration with leading automation manufacturers worldwide, including Microsoft.

OPC provides inter-operability between system components by creating and maintaining open standard specifications. The benefit of following standard specifications in system implementations is greater independence for hardware and software components. This leads to a higher flexibility, better quality, and an overall lower maintenance costs for the system solution. The first standard developed was the Data Access (DA) specification, which therefore has the broadest client support.

The UNICORN OPC Server supports the following three areas:

- The UNICORN OPC DA that gives access to all process data (e.g., real-time values, valve status, process step information, and commands). UNICORN Alarm and Events (AE) Server informs an OPC UA Client application that a system parameter has exceeded an upper or lower limit value.
- The UNICORN Historical DA allows any OPC UA Client application to access the entire batch results generated by UNICORN.
- The UNICORN OPC security controls client access to the UNICORN OPC DA, Alarms and Events, and Historical DA to protect sensitive information and to guard against unauthorized modification of process parameters. This is an important security feature.

1.2 OPC UA standards and versions supported by UNICORN OPC Server

The UNICORN OPC Server supports the following OPC UA standards:

- Data Access and Alarms and Events Server (DA and AE server)
- Historical Data Access (HDA)
- Security

The UNICORN OPC Server supports `Basic256Sha256` and `Aes128_Sha256_RsaOaep` security policies.

The UNICORN OPC Server is certified by the OPC foundation. The profile *Standard 2017 UA Server Profile* was selected and certified, refer to the profile details in the link below:

<https://profiles.opcfoundation.org/profile/1663>

This document requires basic knowledge of UNICORN and the OPC UA Client.

2 UNICORN OPC Server general settings

About this chapter

This chapter provides the required system settings for the UNICORN OPC Server.



IMPORTANT

An OPC UA Client does not need to be installed before performing the steps in this chapter. In this document, UaExpert™ is used as an OPC UA Client for performing the steps. The screenshots are also taken using the UaExpert and defined with a particular system. We recommend uninstalling any previous version of the UNICORN OPC Server before installing the latest version of UNICORN OPC Server 1.3.



IMPORTANT

For a new ÄKTA process™ system, the user must be connected and in control of the system before it is possible to configure the UNICORN OPC Server. For all other system types, the user is not required to be connected to the system.

Installation prerequisites

Install Microsoft .Net Windows Desktop Runtime 6.0.27 x86 and Microsoft ASP.NET Core Runtime 6.0.27 x86 before running the UNICORN OPC Server installer. Click the link below to find the .Net installers:

<https://dotnet.microsoft.com/en-us/download/dotnet/6.0>

UNICORN OPC Server software requirements

Attribute	Specification
Logical processors	Minimum 4 logical processors are recommended
OS version	Windows 10 version 1607 or later (x64) Professional/Enterprise is required
OS language for logged on user	English (U.S.), LCID 1033 is recommended

Attribute	Specification
OS language for system user	English (U.S.), LCID 1033 is recommended
File system format	NTFS is required
Minimum free disk space	Minimum 60 GB
Minimum available RAM	Minimum 8 GB
Page file settings	Minimum 4 GB

In this chapter

Section	See page	
2.1	Install UNICORN OPC Server	9
2.2	UNICORN OPC Server license	11
2.3	Installation log files and server configuration files	13
2.4	Configure UNICORN OPC Server	14
2.5	Configure OPC UA Client and connect to UNICORN OPC Server	18
2.6	Trust certificate	25
2.7	Execute method/manual instruction and raise and acknowledge alarm with the OPC UA Client	28

2.1 Install UNICORN OPC Server

Follow the instructions below to install the UNICORN OPC Server.

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

- | | |
|---|--|
| 1 | Run the UNICORN OPC Server Setup 1.3.exe installation file. |
|---|--|

Note:

If a window appears for installing Microsoft .Net Windows Desktop Runtime 6.0.27 x86 or Microsoft ASP.NET Core Runtime 6.0.27 x86, see [Installation prerequisites, on page 7](#).

- | | |
|---|---|
| 2 | Click Next in the UNICORN OPC Server 1.3 - InstallShield Wizard . |
|---|---|

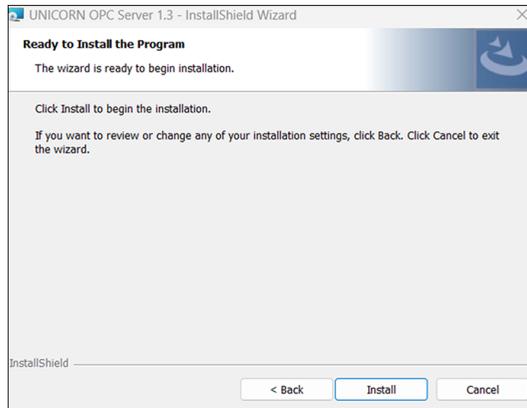


- | | |
|---|---|
| 3 | Agree to the terms in the license agreement, then click Next . |
|---|---|

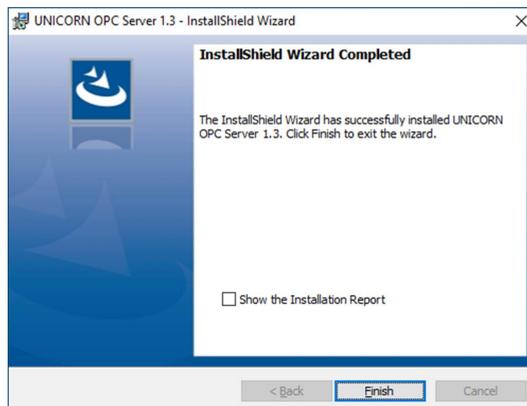


Step Action

- 4 Click **Install** to begin the installation.



- 5 Click **Finish** to close the **InstallShield Wizard**.



2.2 UNICORN OPC Server license

Introduction

It is assumed that the UNICORN OPC Server is installed, and that the UNICORN OPC service is running.

UNICORN OPC Server license (Concurrent) should be generated, and configured using the **Configure e-License** tool.

The license server for UNICORN OPC Server needs to be provided.



IMPORTANT

For the Concurrent license, provide the concurrent license server Hostname or IP address.

The **UNICORN OPC Server.dll.config** file is located at C:\Program Files (x86)\Cytiva\UNICORN OPC Server 1.3\Bin.

Change service startup type to Automatic (Delay Start)

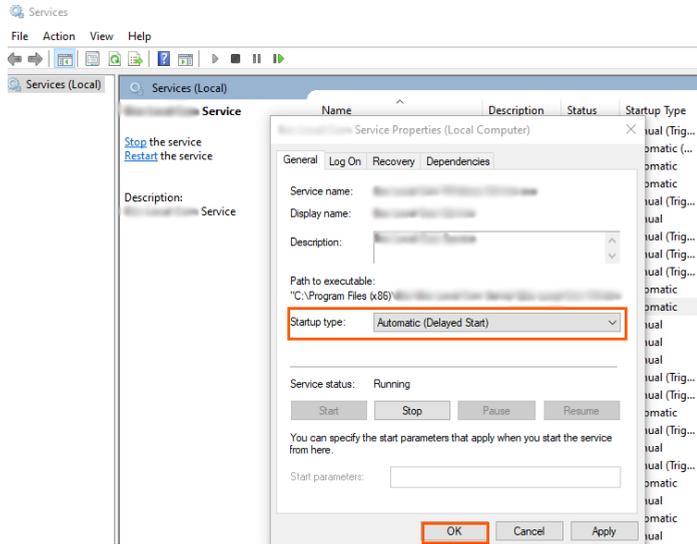
We recommend changing the **Startup Type** of the **UNICORN OPC Server** service to **Automatic (Delay Start)** in the **Services** app in Windows.

Follow the instructions below to perform the change.

Step	Action
1	Open the Services app in Windows.
2	Right-click the UNICORN OPC Server service and click Properties .
3	Change the Startup type to Automatic (Delayed Start) .

Step Action

4 Click **OK**.



2.3 Installation log files and server configuration files

The following table describes the locations of the installation log files and the server configuration files:

Files	Location
Installation log files	C:\ProgramData\Cytiva\UNICORN OPC Server\Logs
UAHDAServer.Config.xml (for HDA Server) UADAAndAEServer.Config.xml (For DA and AE Server) UNICORN OPC Server.dll	C:\Program Files (x86)\Cytiva\UNICORN OPC Server 1.3\Bin

In order to enable more logs, change the value of the tags below in the config file (UNICORN OPC Server.dll).

- `<GatewayLoggerDefaultLogLevel="">`
- `<logger name="GatewayRunLogger">`
- `<logger name="GatewayCoreLogger">`
- `<logger name="WatchdogLogger">`
- `<logger name="DAAndAEAdapterLogger">`
- `<logger name="HDAAdapterLogger">`

In the example below, more logs can be enabled by changing the level value = **Trace**, **Debug**, **Information**, **Warning**, **Error**, or **Critical**.

```
<logger name="GatewayRunLogger">
  <level value="ERROR"/>
  <appender-ref ref="GatewayRunLog"/>
</logger>
```

After changing the level value of any of the tags above, save the config file, and restart UNICORN OPC Server to implement the changes.

Note: When the level value is changed to = **Trace**, **Debug**, **Information**, **Warning**, **Error**, or **Critical**, the size of the log file increases accordingly.

2.4 Configure UNICORN OPC Server

Introduction

The **UNICORN OPC Server Configuration Tool** is used to update the following:

- Database
- License server host name
- Ports
- UNICORN OPC Server status
- Security policies
- Log Levels

Type **UNICORN OPC Server Configuration tool** in **Windows Search** and select **Run as Administrator**.

Use the **UNICORN OPC Server Configuration Tool** window to configure the Database, the License Server host name, the Ports, and also to enable or disable some security policies as detailed in the following sections.

Database Configuration

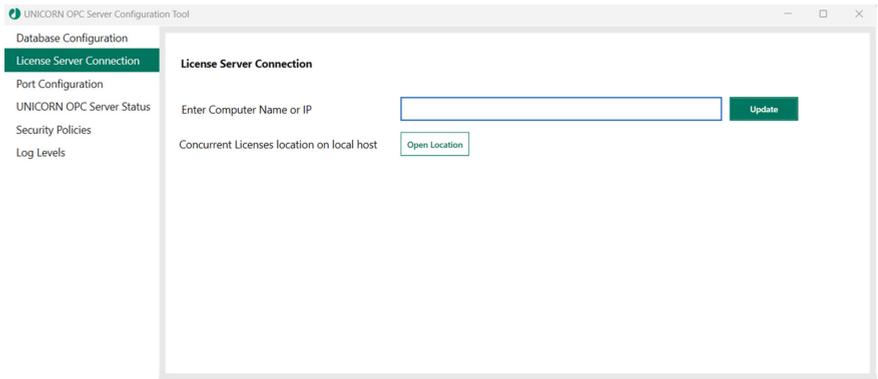
Enter the UNICORN **Database Host Name** or **IP** address and click **Update**. Then the host name or the IP address is added to the configuration file.

License Server Connection

Enter the OPC License Server **Computer Name** or **IP** address and click **Update**. Then the computer name or the IP address is added to the configuration file.

Note: *The License Server for UNICORN OPC Server needs to be provided.*

*For a Concurrent license, the license host value is the concurrent License Server IP address or **Computer name**.*

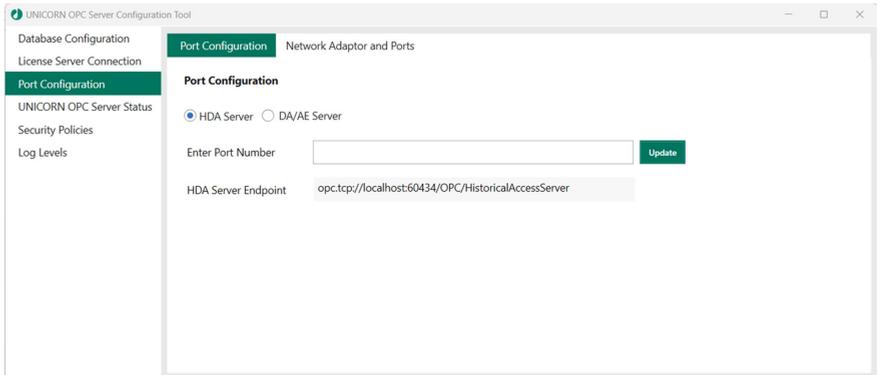


Port Configuration

Port Configuration

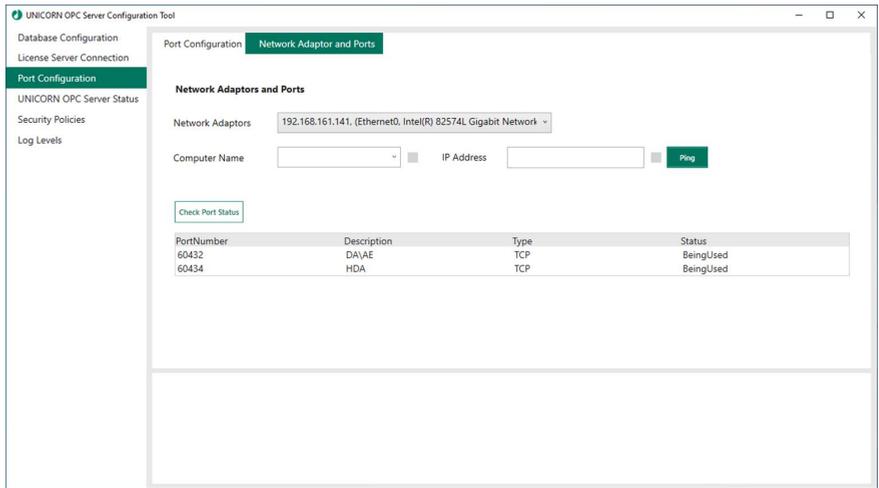
The default Port for the **DA/AE Server** is 60432, and the default Port for the **HDA Server** is 60434.

To change the Port, select **HDA Server** or **DA/AE Server**, type the Port number in the **Enter Port Number** text field, and click **Update**. Then the Port number is added to the configuration file.



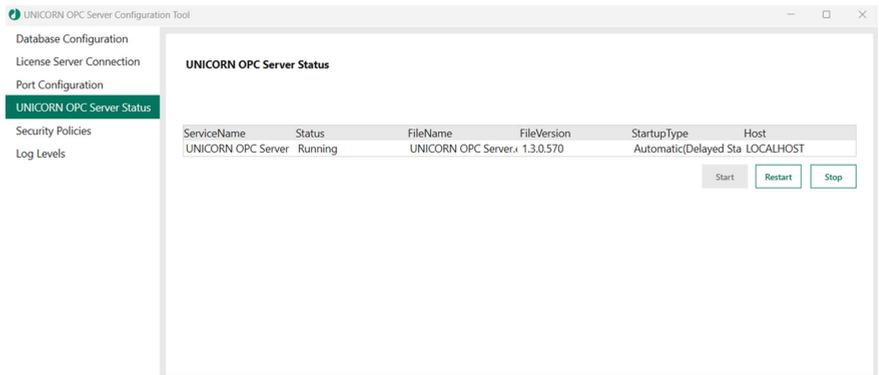
Network Adaptor and Ports

The user can check the **Status/Type** of the **PortNumber** of DA/AE/HDA server, which are assigned in the **Port Configuration** tab on the selected computer name.



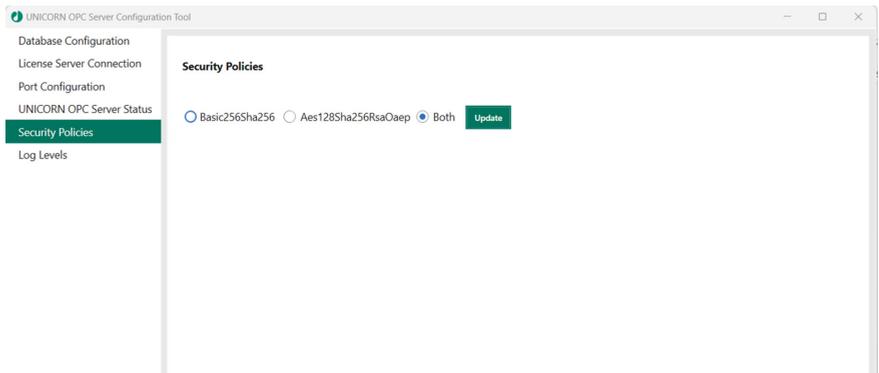
UNICORN OPC Server Status

Use this setting to **Start**, **Restart**, or **Stop** the **UNICORN OPC Server**.



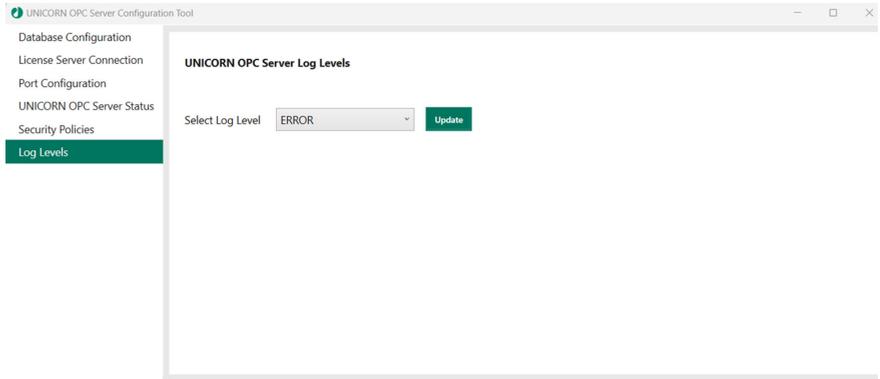
Security Policies

Select any one of the **Security Policies** or **Both** and click **Update**.



Log Levels

Click the drop-down box and **Select Log Level** (ERROR, INFO, WARNING, DEBUG, ALL) to get detailed UNICORN OPC Server logs, then click **Update**.



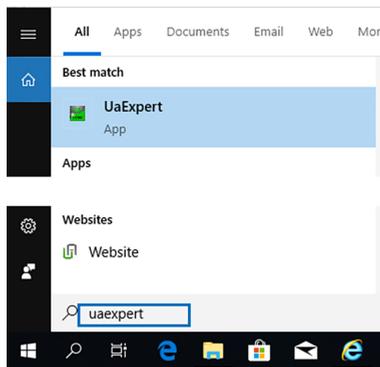
2.5 Configure OPC UA Client and connect to UNICORN OPC Server

Note: In this section, the UaExpert is used as the OPC UA Client for performing the different steps. The screenshots show the UaExpert interface.

Follow the instructions below to configure and connect the OPC UA Client to the UNICORN OPC Server.

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

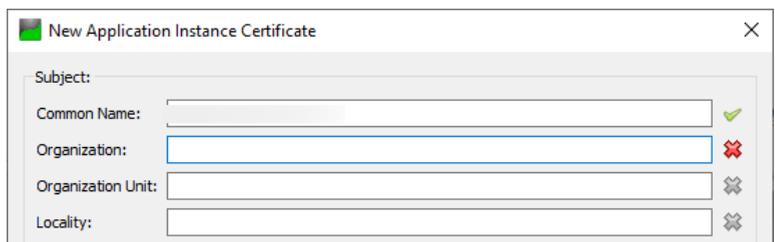
1	Enter <code>uaexpert</code> in the Windows search box and press Enter .
---	--



2	If the following dialog box appears, click OK .
---	--



3	Enter the details for the below fields in the New Application Instance Certificate window.
---	---



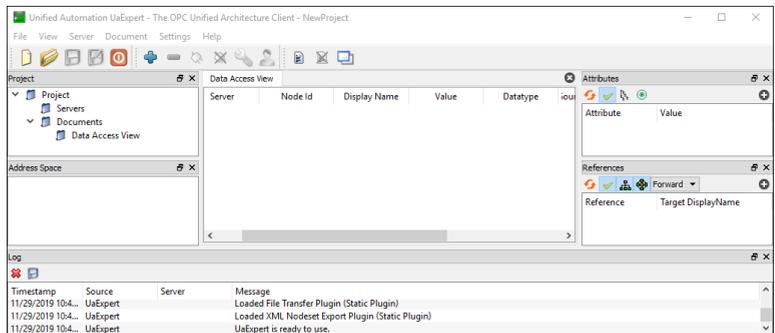
Step Action

4 Click **OK**.



Result:

The following window appears:



Tip:

To avoid the **BadTimeout** error during the connection of the UNICORN OPC Server, or during the browsing and execution of the method call, the user is suggested to increase the **Timeout** value in the OPC UA Client.

The user needs to increase the value of the **General.WatchdogTimeout** if the parameter was subscribed to for monitoring, but the client does not receive the monitoring response (i.e., grayed out for a moment) within the specified **WatchdogTimeout** value.

For reference, the **Timeout** value can be set in the **OPC UAExpert Client**, as below:

Go to **Setting** → **Configure UaExpert**, then set a larger value for the **General.Browse.Timeout** parameter, as seen in the below screenshot:

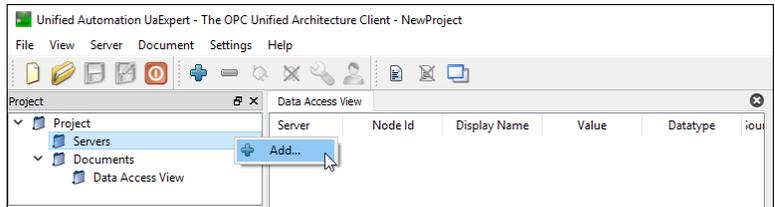
Step Action

Parameter	Value
General.BrowseTimeout	200000
General.CallTimeout	10000
General.ConnectTimeout	10000
General.DiscoveryTimeout	10000
General.InternalServiceCallTimeout	5000
General.PublishTimeout	60000
General.SessionTimeout	3.6e+06
General.WatchdogTimeout	5000
Stack.ThreadPool_Timeout	4294967295

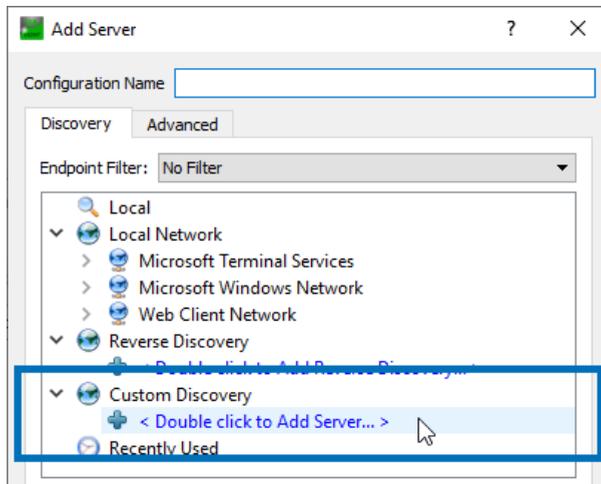
Note:

The **Timeout** value can vary for different OPC UA Clients, and it also depends on the infrastructure.

- 5 Right-click on **Servers** and then click **Add**.



- 6 Double-click on the item as illustrated below:

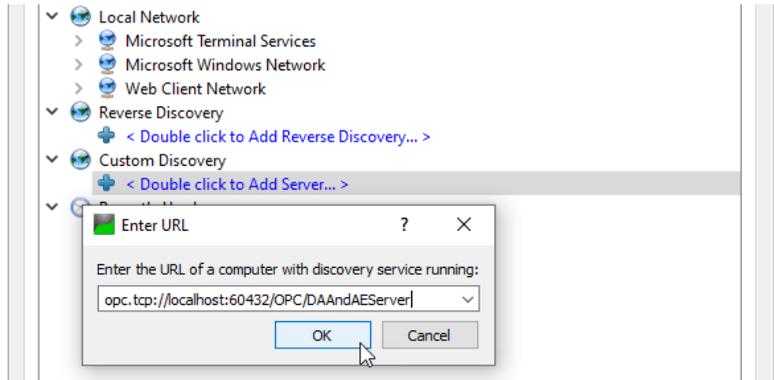


Step	Action
------	--------

7	Enter the IP address/URL of the DA and AE Server/HDA Server and click OK . Below is an example URL of a DA and AE Server.
---	--

Note:

The UNICORN OPC Server URL/Address is available in the UNICORN OPC Server configuration tool.



- Definition of the DA and AE server URL:

opc.tcp://<IP Address/Hostname of the UNICORN OPC Server>:<Port Number of the DAandAE Server>/OPC/<DAandAEServer>

- Definition of the HDA server URL:

opc.tcp://<IP Address/Hostname of the UNICORN OPC Server>:<Port Number of the HDA Server>/OPC/<HistoricalAccessServer>

Example of URL where the OPC Server and the client are on the same machine and default ports are used:

opc.tcp://localhost:60432/OPC/DAAndAEServer

When connecting from a remote client, provide the IP address of the system where the OPC Server is running.

Use an updated port in the URL string if the port has been updated using the OPC Configuration Tool.

Note:

The URL string to add a DA/AE Server or an HDA Server to the OPC UA Clients can be obtained from the configuration files located in *C:\Program Files (x86)\Cytiva\UNICORN OPC Server 1.3\Bin*.

File names:

- DA/AE Server: ***UADAAndAEServer.Config***

Step Action

```

UAADAAndAEServer.Config.xml
74 </TransportQuotas>
75 <ServerConfiguration>
76 <BaseAddresses>
77 <!--<ua:String>https://localhost:60431/OPC/DAAndAEServer</ua:String-->
78 <ua:String>opc.tcp://localhost:60432/OPC/DAAndAEServer</ua:String>
79 </BaseAddresses>
80 <SecurityPolicies>
    
```

- **HDA Server: UAHDAServer.Config**

```

UAHDAServer.Config.xml
73 <SecurityTokenLifetime>3600000</SecurityTokenLifetime>
74 </TransportQuotas>
75 <ServerConfiguration>
76 <BaseAddresses>
77 <!--<ua:String>https://localhost:60433/OPC/HistoricalAccessServer</ua:String-->
78 <ua:String>opc.tcp://localhost:60434/OPC/HistoricalAccessServer</ua:String>
79 </BaseAddresses>
    
```

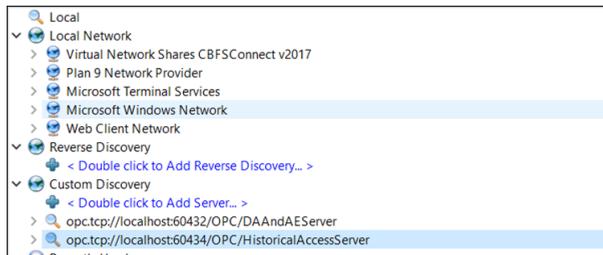
We do not recommend updating the configuration files manually.

- 8 Click **Yes**.



Result:

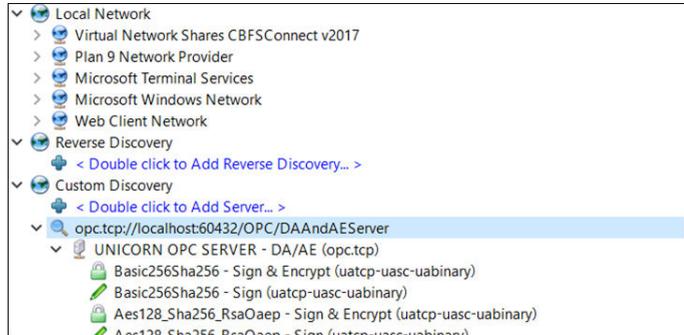
The IP address/URL is now listed under **Custom Discovery**.



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

9	Connect UA Expert with UNICORN OPC Server
---	--

Expand the IP/URL and double-click the highlighted (example) server as illustrated below:



10	Enter UNICORN Username and Password . The users' logons from the OPC UA Client should be part of only one access group.
----	---

Note:

We do not recommend disabling the password in UNICORN while using UNICORN OPC Server.

If the user disables the password in UNICORN, the user should provide a password identical to their username when logging on to the OPC UA client.

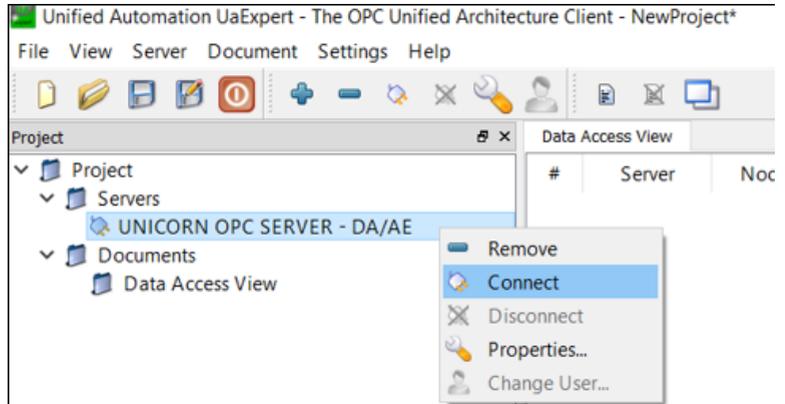
Note:

We do not recommend changing the logged on users from OPC UA client without closing the session completely.

Example: If the default user has logged on to the OPC UA client, do not switch to another user until the session is completely disconnected

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

- | | |
|----|--|
| 11 | Go back to the Unified Automation UaExpert window and right-click on OPC DA and AE server and the click Connect . |
|----|--|



Note:

The connected OPC UA user can be identified by **OPC_Hostname** displayed in system control, in UNICORN, and system logs.

2.6 Trust certificate

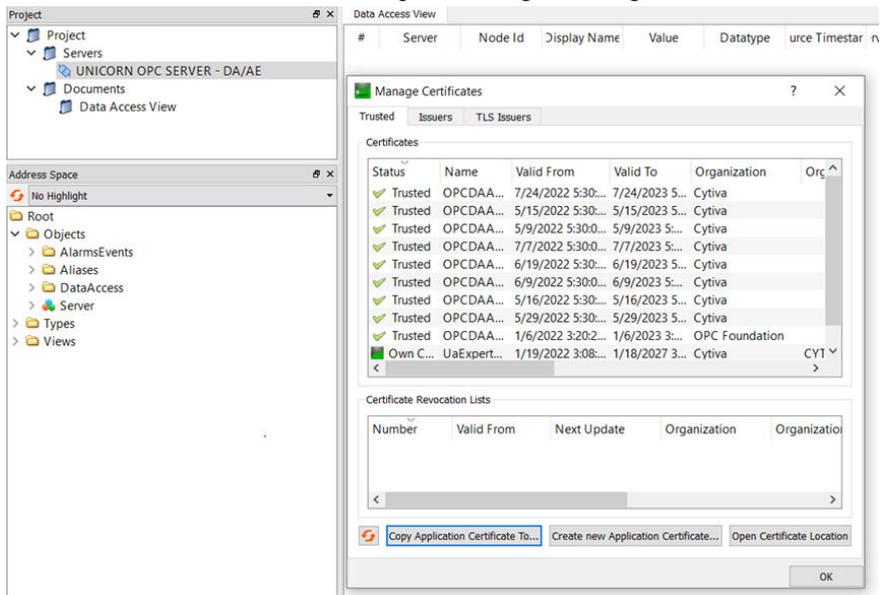
An initial attempt to securely connect to an UNICORN OPC Server from the Workflow server can fail due to a missing client certificate with the following error message:

Bad Security Check Failed.

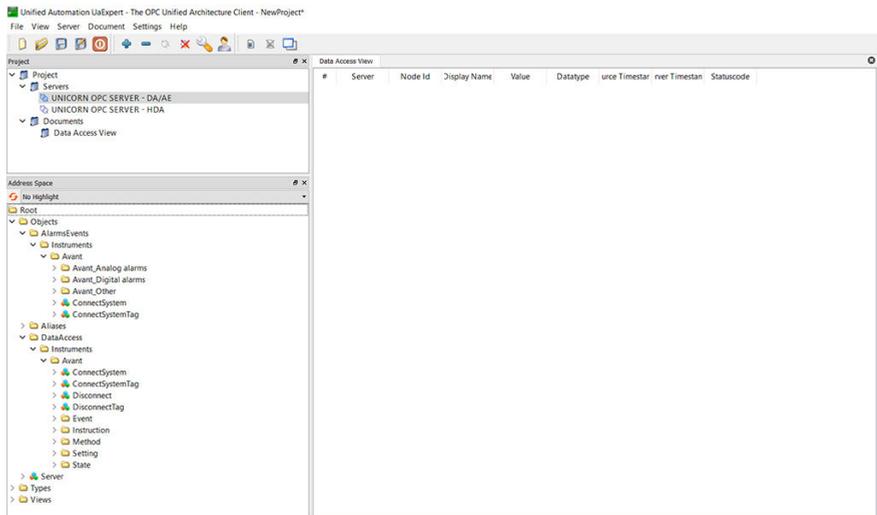
Once a certificate is created, you can establish a connection to the UNICORN OPC Server. The process is described below:

1. The certificate is sent to the server as a part of the connection request.
2. The UNICORN OPC Server returns the server certificate during the connection exchange.
3. The UNICORN OPC Server does not trust the client certificate initially and places the certificate in `C:\ProgramData\Cytiva\UNICORN OPC Server\pki\rejected\certs`.
4. Cut the certificate from the above location and paste it to the location below:
`C:\ProgramData\Cytiva\UNICORN OPC Server\pki\trusted\certs`.

To check the details of the certificates, go to **Settings → Manage Certificates**.



The address space appears after connection to the server.



UNICORN OPC Server creates a self-signed certificate by default.

Note:

- *When UNICORN OPC Server certificate has expired during an ongoing run then there will not be a communication break down from the current session.*

An ongoing run will not be impacted if UNICORN OPC Server certificate has.

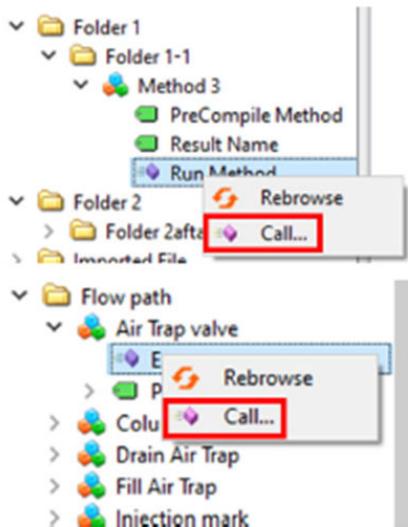
- *Certificate renew should be done when no connection to UNICORN OPC Server exists.*
- *When a user creates a new connection we recommend that the user manually deletes both OPCDAAndAEServer and HDAServer certificates from the path below and restart the UNICORN OPC Server so that OPC stack validates and creates a new certificate:*

```
C:\ProgramData\Cytiva\UNICORN OPC Server\pki\own\certs
```

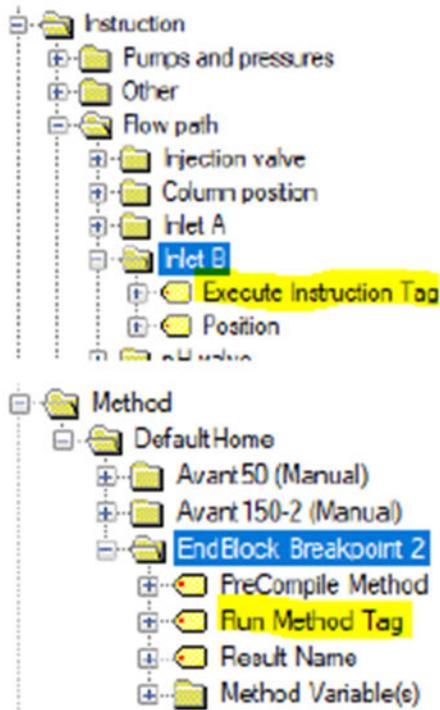
```
C:\ProgramData\Cytiva\UNICORN OPC Server\pki\trusted\certs
```

Production Client Settings

For the OPC UA Client that supports **Method Call / Method services**, the UNICORN method is executed by clicking **Call...**, as shown in the screenshots below.



For the OPC UA Client that does not support **Method Call / Method services**, the UNICORN method cannot be executed by clicking **Call...**. Instead, **Execute Instruction Tag**, **Execute Setting Tag**, or **Run Method Tag** can be clicked to execute the UNICORN method, as shown in the example screenshots below.



2.7 Execute method/manual instruction and raise and acknowledge alarm with the OPC UA Client

Introduction

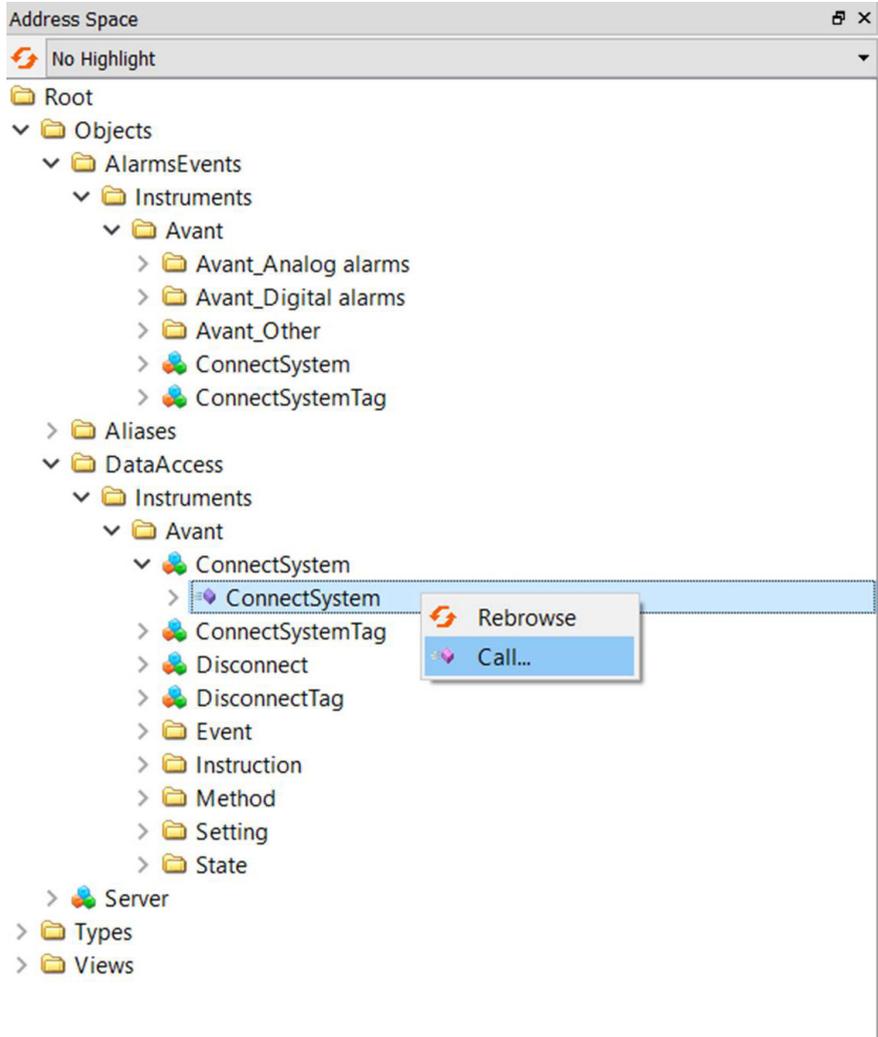
This section provides information regarding how to use the OPC UA Client to execute a method or a manual instruction, and how to raise and acknowledge an alarm.

In this section

Section	See page
2.7.1 Connect OPC Client with Instrument	29
2.7.2 Execute method	30
2.7.3 Execute manual instruction	32
2.7.4 Raise and acknowledge an alarm	34
2.7.5 Disconnect OPC Client with Instrument	38

2.7.1 Connect OPC Client with Instrument

Call the **ConnectSystem** method to connect the OPC Client to the Instrument.

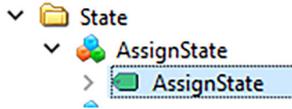


2.7.2 Execute method

Follow the instructions below to execute a method using the OPC UA Client.

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

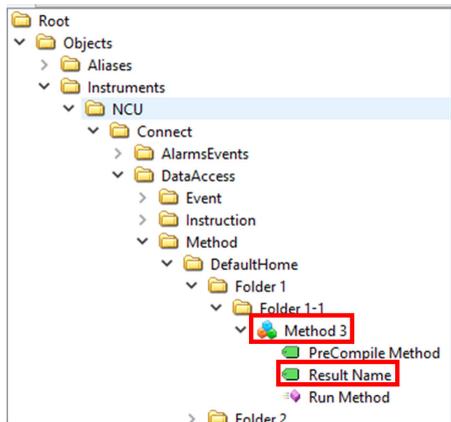
- | | |
|---|---|
| 1 | Subscribe to the AssignState tag, and set the value to 1 to connect the OPC in control mode. |
|---|---|



Note:

A value of 0 means the OPC is connected in view mode.

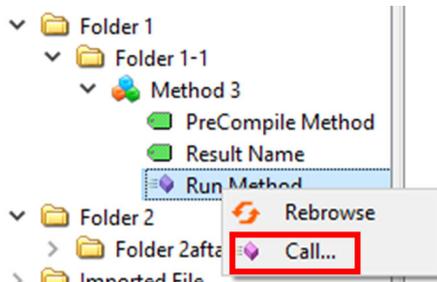
- | | |
|---|---|
| 2 | Navigate to the Method , then drag and drop Method 3 , or the Result Name tag. |
|---|---|



- | | |
|---|---|
| 3 | Enter the Result Name for the file, for example, OPCTestRun . |
|---|---|

#	Server	Node Id	Display Name	Value
1	UNICORN OPC SERVER - DA/AE	NS3\$String1[NCU]AssignState:DefaultAssignState	AssignState	1
2	UNICORN OPC SERVER - DA/AE	NS3\$String1[NCU]Method 3:Folder 1-1:PreCompile Method	PreCompile Method	
3	UNICORN OPC SERVER - DA/AE	NS3\$String1[NCU]Method 3:Folder 1-1:Result Name	Result Name	"1.0?"; <Method Name="Method 3" Id="777"; <Status>Succeed</Status></...>

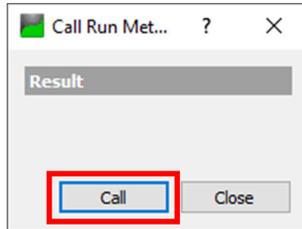
- | | |
|---|---|
| 4 | Right-click Run Method , then click Call... |
|---|---|



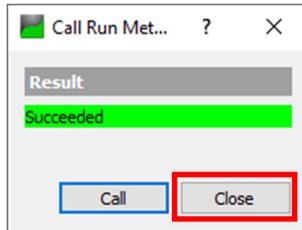
Step	Action
------	--------

Result:

Click **Call...** in the **Call Run Met...** window that appears.

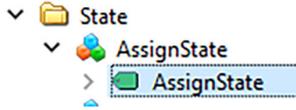
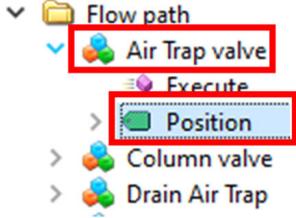
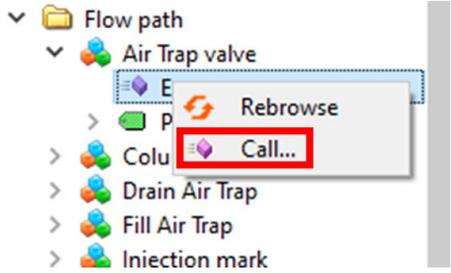
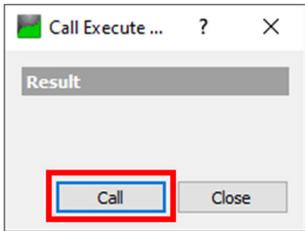


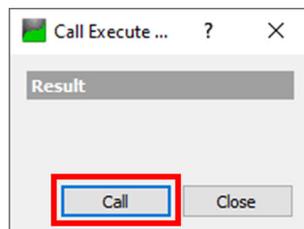
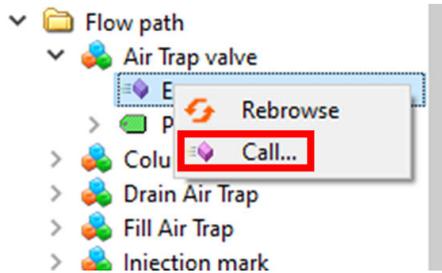
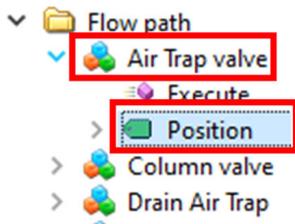
- 5 If the call is successful when **Succeeded** appears, then click **Close**.



2.7.3 Execute manual instruction

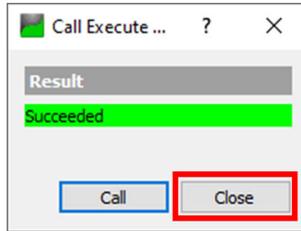
Follow the instructions below to execute a manual instruction using the OPC UA Client.

Step	Action
1	<p>Subscribe to the AssignState tag, and set the value to 1 to connect the OPC in control mode.</p> 
	<p>Note: A value of 0 means the OPC is connected in view mode.</p>
2	<p>Subscribe to a tag, for example, AirTrap Valve, then drag and drop the Position tag. Set the value to 0 for bypass or 1 for inline.</p> 
3	<p>Right-click Execute, then click Call...</p> 
	<p>Result: Click Call... in the Call Execute ... window that appears.</p> 



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

4	If the call is successful when Succeeded appears, then click Close .
---	--



Step Action

- 3 Select and right-click the alarm, then click on **Acknowledge** to acknowledge the alarm. The acknowledged alarm will now be logged under the **Events** tab.



Note:

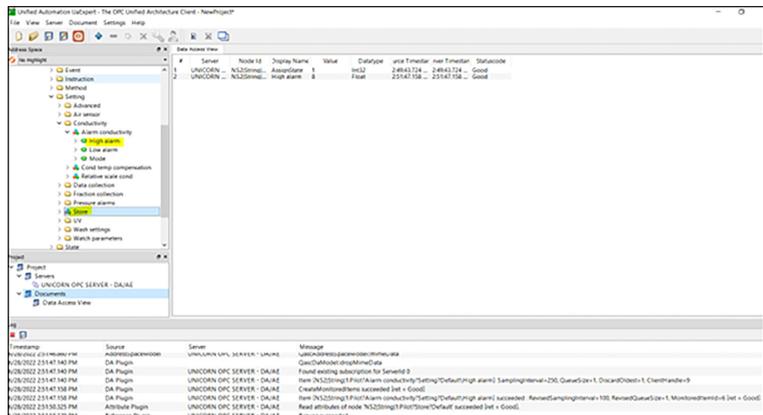
The alarms disappear from the queue once they are acknowledged.

Raise an analog alarm for **Alarm conductivity tag**

Follow the instructions below to raise an analog alarm for the **Alarm conductivity** tag using the OPC UA Client.

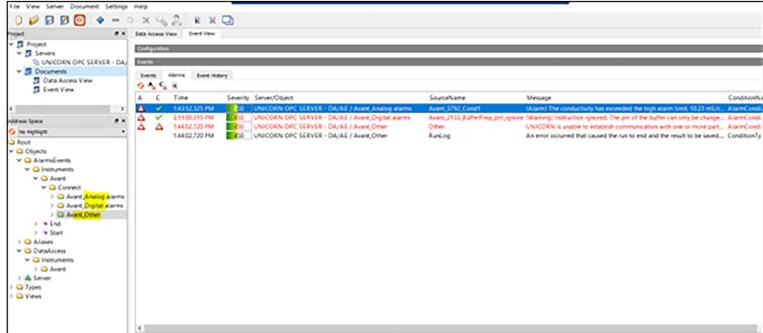
Step Action

- 1 In the /State/AssignState file path under the **DA** server, add the **AssignState** tag. In the /Setting/Alarm conductivity file path, add the **High alarm** and **Execute** tags.

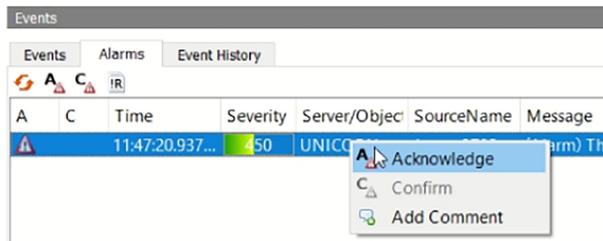


Step Action

- 2 Click the **Alarms** tab to see the generated analog alarm.



- 3 Select and right-click the alarm, then click on **Acknowledge** to acknowledge the alarm. The acknowledged alarm will now be logged under the **Events** tab.



Note:

The alarms disappear from the queue once they are acknowledged.

Raise a digital alarm for BufferPro pH tag

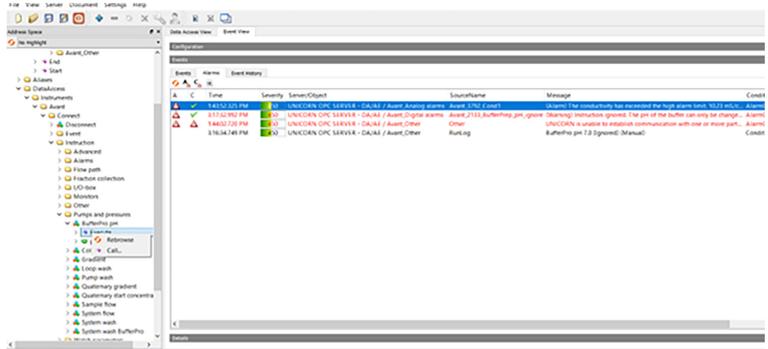
Follow the instructions below to raise an digital alarm for the **BufferPro pH** tag using the OPC UA Client.

Step Action

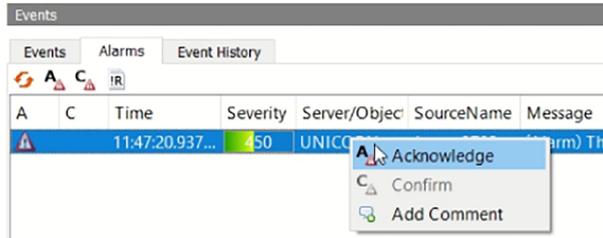
- 1 In the /State/AssignState file path under the **DA** server, add the **AssignState** tag. In the /Instruction/Pumps and pressures file path, add the **BufferPro pH** and **Execute** tags.

Step Action

- 2 Click the **Alarms** tab to see the generated digital alarm.



- 3 Select and right-click the alarm, then click on **Acknowledge** to acknowledge the alarm. The acknowledged alarm will now be logged under the **Events** tab.



Note:

The alarms disappear from the queue once they are acknowledged.

2.7.5 Disconnect OPC Client with Instrument

Subscribe to the **Disconnect** tag, and set the value to 1 to disconnect the OPC from the Instrument.

The **Disconnect** method call will also disconnect the OPC from the instrument.

The screenshot shows the Unified Automation UaExpert interface. The 'Data Access View' window displays the following data:

#	Server	Node Id	Display Name	Value	Datatype	urce Timestar	ver Timestar	Statuscode
1	UNICORN ...	NS2[String]...	Disconnect	1	Int32	5:56:46.644 ...	5:56:46.644 ...	GoodNoData

The 'Project' tree on the left shows the following structure:

- Project
 - Servers
 - UNICORN OPC SERVER - DA/AE
 - UNICORN OPC SERVER - HDA
 - Documents
 - Data Access View
- Address Space
 - No Highlight
 - Root
 - Objects
 - AlarmsEvents
 - Instruments
 - Avant
 - Avant_Analog alarms
 - Avant_Digital alarms
 - Avant_Other
 - ConnectSystem
 - ConnectSystemTag
 - Aliases
 - DataAccess
 - Instruments
 - Avant
 - ConnectSystem
 - ConnectSystemTag
 - Disconnect
 - DisconnectTag
 - Event
 - Instruction
 - Method

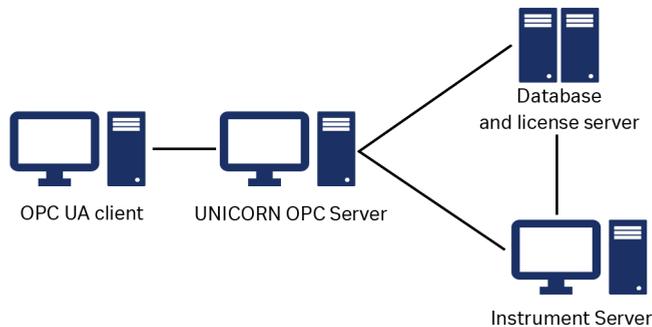
3 Deployments

Introduction

This chapter describes the various deployments by which the UNICORN OPC Server can be connected to the OPC UA Client.

Distributed environment (single system)

The following image shows a single system deployment.



Distributed environment (multisystem)

The UNICORN OPC Server has passed the 72 hour run, required by the OPC foundation. We recommend using a maximum of 5 users or a maximum of 5 systems if used in a distributed/multisystem environment with a monitoring load of 5K tags.



IMPORTANT

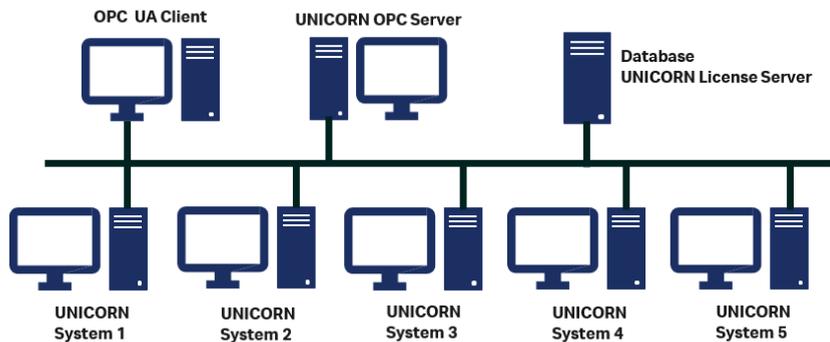
We do not recommend installing the UNICORN OPC Server on ReadyToProcess™ 25.

Note: *The user can connect more than five systems with one UNICORN OPC Server, but performance or connectivity issues could be encountered.*

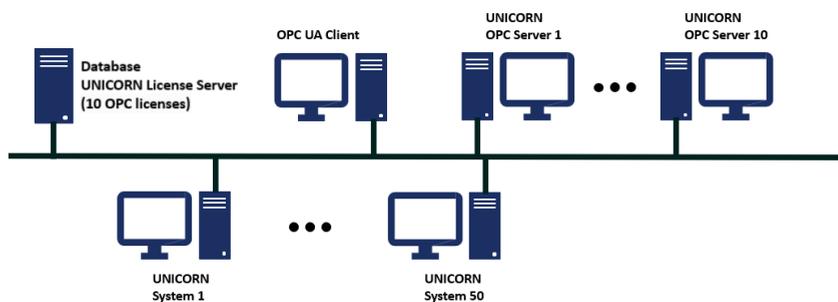
Note: *Do not connect multiple OPC UA Clients (from the same or different computer) to the same system. If multiple users or the same user needs to connect to same system, then log off the current OPC UA user and log on with the user that needs to view the data. Disconnecting one user will disconnect all the other users who are connected to the same system.*

Note: *If the same user wants to connect to different instruments from two different computers using the UNICORN OPC Server, we recommend the following scenarios:*

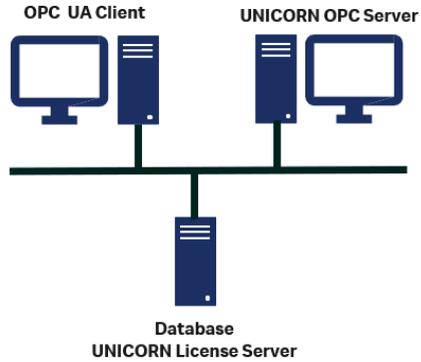
- *The same and unique user id should be used when connecting to a specific instrument, so one dedicated user should be used for each instrument.*
- *The same and unique user id should be used when connecting from a specific client.*
- *Alternatively, make sure all the connections are performed from the same client.*
- *Use one UNICORN OPC Server for each instrument, but this can increase the cost of a license.*



The below image shows the recommended setup for the distributed environment with more than five systems.



The below image shows the Historical Data Access (HDA) Server connection. In this HDA Server connection, the user can access the UNICORN OPC HDA address space, audit trail, HDA XML format definition, and so on, without connecting to the instrument.



Note: *The UNICORN Database and license server can be installed on a different machine.*

4 UNICORN OPC Data Access address space

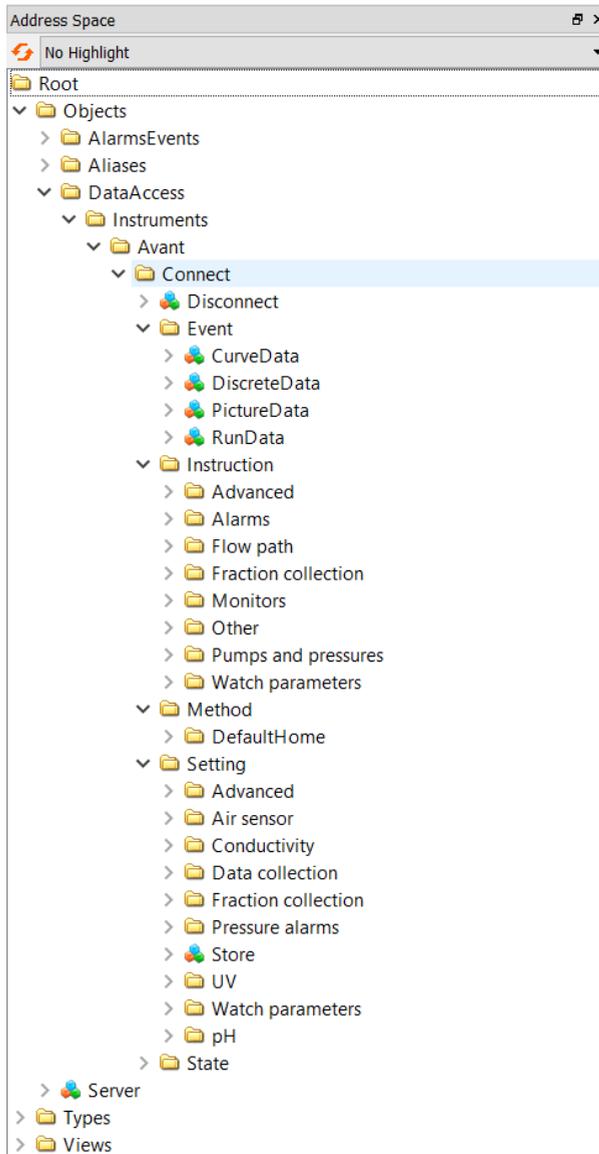
About this chapter

This chapter provides information about the UNICORN OPC Historical Data Access (HDA) address space, method execution, settings, run data, picture data, and so on.

The OPC DA address space is a real-time representation of the process control data for the connected instruments.

Data Access address space view

The following screenshot shows the DA address space view.



Tip: *The address space and the tags depend on the instrument configurations, connected systems, and the user logged on. The OPC server tags contain a username that is included in the Node Id (tag path). The tags generated under the **Event**, **Instruction**, **Setting**, and **State** folders include the user that is logged on. For example, the default user is logged on into the OPC UA*

Client in all the screenshots in the sections of this chapter. See the screenshots in the sections of this chapter to see how the tags are built. The tags are configured by a user during the initialization of the production client. The same user needs to log into the production client to access the configured tags.

In this chapter

Section		See page
4.1	Run data and Picture data	45
4.2	Trend data	48
4.3	Manual instructions	51
4.4	Method execution	52
4.5	State	67
4.6	Recommendations	79

4.1 Run data and Picture data

Introduction

UNICORN run data are available via the `EVENT\Rundata` branch, and can be analog or digital depending on the data type. Picture data are available via `EVENT\PictureData`.

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
Event\Rundata\Accumulated time
```

```
Event\Rundata\System flow
```

```
Event\Rundata\BufferPro pH
```

```
Event\Analog Picture\UV1
```

```
Event\Digital Picture\Frac Comp
```

Properties for Run data

Properties for Run data are: (Event\Rundata\System flow)

The screenshot shows a window titled 'Attributes' with a toolbar containing icons for refresh, check, help, and a plus sign. The main area is a table with two columns: 'Attribute' and 'Value'. The 'NodeId' attribute is expanded to show a list of sub-attributes. Below the table is a 'References' section with a toolbar and a table of references.

Attribute	Value
▼ NodeId	ns=2;s=1:AvantSystem?RunData?Default\System flow
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?RunData?Default\System flow
NodeClass	Variable
BrowseName	2, "System flow"
DisplayName	"" , "System flow"
Description	"" , "System flow"
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
▼ Value	
SourceTimestamp	03-03-2022 12:25:35.725
SourcePicoseconds	0
ServerTimestamp	03-03-2022 12:25:35.724
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	0
▼ DataType	Float
NamespaceIndex	0
IdentifierType	Numeric
Identifier	10 [Float]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EURange
HasProperty	EngineeringUnits
HasTypeDefiniti...	AnalogItemType

Properties for Picture data

Properties for Picture data are: (Event\Analog Picture\UV1)

The screenshot shows a window titled 'Attributes' with a toolbar at the top containing icons for refresh, check, help, and a green circle. The main area is a table with two columns: 'Attribute' and 'Value'.

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?PictureData?Default\UV 1
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?PictureData?Default\UV 1
NodeClass	Variable
BrowseName	2, "UV 1"
DisplayName	"" , "UV 1"
Description	"" , "UV 1"
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 12:29:11.773
SourcePicoseconds	0
ServerTimestamp	03-03-2022 12:29:11.772
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	0
Data Type	Float
NamespaceIndex	0
IdentifierType	Numeric
Identifier	10 [Float]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Below the 'Attributes' section is a 'References' section with a toolbar containing icons for refresh, check, help, and a green circle, followed by a dropdown menu set to 'Forward'. The main area is a table with two columns: 'Reference' and 'Target DisplayName'.

Reference	Target DisplayName
HasProperty	EURange
HasProperty	EngineeringUnits
HasTypeDefiniti...	AnalogItem Type

4.2 Trend data

Introduction

The UNICORN trend data are available via the `EVENT\Curves`. These data have the highest resolution. **ANALOGTREND** data are always an array of values.

BINARYTREND data

BINARYTREND contains textual trending data:

Data	Description
CurrentMethod	Name of the current method run.
CurrentResult	Name of the current result to which the method run is saved.
CurrentScouting	Scouting number.
CurrentBlock	Name of the block currently executing.
RunLog	Messages from the current run.
Fractions	Current fraction mark.
Injections	Current injection mark.

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
Event\CurveData\UV1
```

```
Event\DiscreteData\Injections
```

Properties for Analog data

Properties for Analog data are: (Event\CurveData\UV1)

Attribute	Value
NodeId	ns=2;s=1:Avant 23?CurveData?Default\UV 1
NamespaceIndex	2
IdentifierType	String
Identifier	1:Avant 23?CurveData?Default\UV 1
NodeClass	Variable
BrowseName	2, "UV 1"
DisplayName	"", "UV 1"
Description	"", "UV 1"
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	1/25/2024 5:59:33.091 PM
SourcePicoseconds	0
ServerTimestamp	1/25/2024 5:59:33.087 PM
ServerPicoseconds	0
StatusCode	GoodNoData (0x00a50000)
Value	Float Array[1]
[0]	0
DataType	Float
NamespaceIndex	0
IdentifierType	Numeric
Identifier	10 [Float]
ValueRank	1 (OneDimension)
ArrayDimensions	UInt32 Array[1]
[0]	0
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	100

Properties for Binary data

Properties for Binary data are: (Event\DiscreteData\Injections)

The screenshot displays two windows from a software interface. The top window, titled 'Attributes', shows a tree view of properties for a binary data point. The bottom window, titled 'References', shows a table of references.

Attributes Window:

Attribute	Value
▼ NodeId	ns=2;s=1:AvantSystem?DiscreteData?Default\Injections
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?DiscreteData?Default\Injections
NodeClass	Variable
BrowseName	2, "Injections"
DisplayName	"" , "Injections"
Description	"" , ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
▼ Value	
SourceTimestamp	03-03-2022 06:28:38.625
SourcePicoseconds	0
ServerTimestamp	03-03-2022 12:34:33.888
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
▼ DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

References Window:

Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DataItem Type

4.3 Manual instructions

Introduction

The UNICORN manual instructions from the **Manual Instructions** box in **System Control** are available in the **INSTRUCTION** branch. To execute a manual instruction, the user can either click on **Call...** to execute a method, or the **Execution Instruction Tag** can be used by providing the parameter value 1 to execute an item, or providing the parameter value 0 to ignore an instruction.

If the instruction in UNICORN uses text, it is possible to get that text from the EU type.

During method runs, manual instruction parameters will be updated with current values, which means that flow changes will be reflected in the flow parameter.

Execute Parameter

Execute parameter is a Method call without any arguments.

Note: *The instructions can only be executed if the module is in control of UNICORN system.*

Quality is good on successful execution of the instruction, which means that the corresponding manual command is successfully submitted.

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\Instruction\Pumps and pressures\System flow\Execute
```

4.4 Method execution

Introduction

The methods are visible in the OPC UA Client after they are created in the UNICORN **Method** module.

A UNICORN method can be run via the **METHOD** branch.

The **METHOD** branch is dynamic. Whenever a new method is created or an existing method is modified, the **METHOD** branch has to be browsed again to see the changes and the **METHOD** branch has to reconfigure the items to use. If a method is deleted, it is removed from the address space.

The tags generated under the **Method** folder do not include the user logged on.

Each method file contains several items.

Note: *Monitoring Method tag values will not update if the user re-browsed the main **METHOD** folder. To get the updated value reflected in the OPC UA Client, the user has to browse the method that they created.*

In this section

Section	See page
4.4.1 Scout Run Index	53
4.4.2 Run Method	55
4.4.3 Result Name	56
4.4.4 Batch ID	58
4.4.5 StartNotes	60
4.4.6 MethodNotes	61
4.4.7 PreCompile	62
4.4.8 Questions	63
4.4.9 Variables	65

4.4.1 Scout Run Index

Introduction

The **Scout Run Index** is used to include any scout runs. When a client writes 0, all scout runs will be included in the method run, unless the excluded flag is set for any particular run. Entering any particular scout run number will include only that run in the method run. The data type is **Int32**.

The object is always **CurrentRead** and **CurrentWrite**. The object is writable only when in control of the system. Executing the method stores the text in the result.

If the user wants to insert a value for the scout variable, then the user needs to first insert a value for the **Scout Run Index** tag to specify the index. When the value is changed for the **Scout Run Index** tag, the user needs to re-browse the method to see the scout variable value for the particular **Scout Run Index**.

Implemented properties

The implemented scout run index properties are as follows:

Attributes	
Attribute	Value
NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\ScoutRunIndex
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\ScoutRunIndex
NodeClass	Variable
BrowseName	2, "ScoutRunIndex"
DisplayName	"" , "ScoutRunIndex"
Description	"" , ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 00:00:00.000
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:09:38.873
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	0
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false
References	
Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DataItem Type

Location (<=> **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem (Manual)
\SampleTestMethod1\Scout Run Index
```

4.4.2 Run Method

The available methods are read from the shared directory of the current user, OPC user, or all users, depending on the setup. **Run Method** is a method call with no arguments.

When a client writes **3** to the state command to end of the method run, the current run state is checked to see if the result is to be saved or not. Other run states imply that it is a manual run being executed, and the result is saved in the [SystemName] (Manual) folder.

Location (<=> **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem (Manual) \SampleTestMethod  
\Run Method
```

4.4.3 Result Name

Introduction

The data type is `String`. Result name objects are both **CurrentRead** and **CurrentWrite**, but only readable from `CACHE`. If a result name is not set in OPC, then UNICORN generates a unique result name when the method starts.

Note: *The methods can only be run if the module is in control of the UNICORN system.*

Implemented properties

The implemented result name properties are as follows:

Attributes	
Attribute	Value
▼ Nodell	ns= 2; s= 1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\Result Name
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\Result Name
NodeClass	Variable
BrowseName	2, "Result Name"
DisplayName	"" "Result Name"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
▼ Value	
SourceTimestamp	03-03-2022 13:21:48.588
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:21:48.586
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
▼ DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false
References	
Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DatalItemType

4 UNICORN OPC Data Access address space

4.4 Method execution

4.4.3 Result Name

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem1 (Manual)
\SampleTestMethod1\Result Name
```

4.4.4 Batch ID

Introduction

The data type is `String`. Batch ID objects are both **CurrentRead** and **CurrentWrite**. If a Batch ID is not set in OPC, then UNICORN generates a unique Batch ID when the method starts.

Note: *The methods can only be run if the module is in control of the UNICORN system.*

Implemented properties

The implemented Batch ID properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\BatchId
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\BatchId
NodeClass	Variable
BrowseName	2, "BatchId"
DisplayName	"" , "BatchId"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 00:00:00.000
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:23:49.955
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DataltemType

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

4 UNICORN OPC Data Access address space

4.4 Method execution

4.4.4 BatchID

```
\Method\DefaultHome\AvantSystem1 (Manual)  
\SampleTestMethod1\Batch ID
```

4.4.5 StartNotes

Introduction

The **StartNotes** is used to read and write the start notes. The data type is `String`. **StartNotes** objects are both **CurrentRead** and **CurrentWrite**. The object is writable only when in control of the system. Executing the method stores the text in the result.

Implemented properties

The implemented start notes properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\StartNotes
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\StartNotes
NodeClass	Variable
BrowseName	2, "StartNotes"
DisplayName	"" "StartNotes"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 13:27:26.587
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:27:26.584
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DataItem Type

The following is an example location ($\langle = \rangle$ **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem1 (Manual)
\SampleTestMethod1\Start Notes
```

4.4.6 MethodNotes

Introduction

The **MethodNotes** is used to read the method notes. The data type is `String`.

MethodNotes objects are **CurrentRead**.

Implemented properties

The implemented method notes properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\Method Notes
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\Method Notes
NodeClass	Variable
BrowseName	2, "Method Notes"
DisplayName	"" , "Method Notes"
Description	"" , ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 00:00:00.000
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:29:02.527
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	TestMethod
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	Definition
HasProperty	ValuePrecision
HasTypeDefiniti...	DataItem Type

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem1 (Manual)
\SampleTestMethod1\Method Notes
```

4.4.7 PreCompile

The data type is always `String`. This forces UNICORN OPC to compile the method and check if it is runnable on the system. If it is runnable on the system, the result will be an xml string containing `Succeed` as the status. If it is not runnable on the system, the string will contain an error text.

Implemented properties

The implemented method compile properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\PreCompile Method
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)\PreCompile Method
NodeClass	Variable
BrowseName	2, "PreCompile Method"
DisplayName	"" "PreCompile Method"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 13:30:42.742
SourcePicoseconds	0
ServerTimestamp	03-03-2022 13:30:42.740
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	<?xml version="1.0"?><Method Name="SampleTestMethod" Id="1172"><Status>Succeed</Status></Method>
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target	DisplayName
HasProperty	Definition	
HasProperty	ValuePrecision	
HasTypeDefiniti...	DataItem Type	

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem (Manual)
\SampleTestMethod1\PreCompile Method
```

4.4.8 Questions

Introduction

The **QUESTIONS** branch is used to read and write the start protocol questions. One object is created for each question in a start protocol. Objects are added to the **QUESTIONS** branch for every method. The canonical data type is either **String** or **Int32**, depending on the question type.

Table 4.1: Question types and data types

Data type	Question type
NoReply	Int32
Entryfield	String
Multiple choice	Int32
Integer	Int32
Float	Float

Integer and **Float** can have minimum and maximum ranges. The writing out of range is not allowed. Empty strings are valid unless the mandatory flag is set. Authorized questions are allowed without supplying any signature. Also, it is possible to run methods without answering all the questions, unless the mandatory flag is set.

Question objects are **CurrentRead** and **CurrentWrite**. The object is writable only when in control of the system. Executing the method stores the answers in the result.

Implemented properties

The implemented properties are as follows:

Attribute	Value
▼ NodeId	ns=2;s=1:AvantSystem?SampleTestMethod?AvantSystem (Manual)?Question(s)\Enter Numeric values
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleTestMethod?AvantSystem (Manual)?Question(s)\Enter Numeric values
NodeClass	Variable
BrowseName	2, "Enter Numeric values"
DisplayName	"" "Enter Numeric values"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
▼ Value	
SourceTimestamp	03-03-2022 13:38:39.003
SourcePicoSeconds	0
ServerTimestamp	03-03-2022 13:41:56.562
ServerPicoSeconds	0
StatusCode	Good (0x00000000)
Value	1
▼ DataType	
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EURange
HasProperty	EngineeringUnits
HasTypeDefiniti...	AnalogItem Type

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\Method\DefaultHome\AvantSystem (Manual)
\SampleTestMethod1\Question(s)\Enter Numeric Values
```

4.4.9 Variables

Introduction

The **VARIABLES** branch is used to read and write the method variables. One object is created for each method variable. Objects are added to the **VARIABLES** branch for every method, if the **StartProtocol** for the method supports the variable value change. The variables on the following are not supported:

- Frac-950 Fractionation instruction parameters
- Frac-950 Peak Fractionation parameters

Variables objects are **CurrentRead** and **CurrentWrite**. The object is writable only when in control of the system. Executing the method uses the variable settings during the method run.

Implemented properties

The implemented start protocol variables general properties are as follows:

Attribute	Value
Node	ns=2;s=1:AvantSystem?SampleMethod?AvantSystem (Manual)?Method Variable(s)\Inlet B
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?SampleMethod?AvantSystem (Manual)?Method Variable(s)\Inlet B
NodeClass	Variable
BrowseName	2, "Inlet B"
DisplayName	"" "Inlet B"
Description	"" ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 15:19:57.754
SourcePicoSeconds	0
ServerTimestamp	03-03-2022 15:19:57.749
ServerPicoSeconds	0
StatusCode	Good (0x00000000)
Value	0
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

4 UNICORN OPC Data Access address space

4.4 Method execution

4.4.9 Variables

```
\Method\DefaultHome\AvantSystem(Manual)  
\SampleTestMethod1\Method Variable(s)\Inlet B
```

4.5 State

Introduction

The **STATE** branch contains important state information for the system.

In this section

Section		See page
4.5.1	AssignState	68
4.5.2	RunState	70
4.5.3	Command	72
4.5.4	InstrumentStatus	74
4.5.5	InstrumentStatusConnection	76
4.5.6	UserInControl	78

4.5.1 AssignState

Introduction

AssignState enables the client to change the assign mode to the UNICORN system. The default assign mode is **View**, which is a mode where the client can monitor the **EVENT** and **STATE** branch items. By changing the mode to **Control**, it is also possible to execute instructions, run methods, and save system settings.

Object values

The defined assign state values are as follows:

Value	State
0	View
1	Control
2	None
3	Disconnected

The quality is good if a correct **AssignState** is written to the object, otherwise it is bad.

Implemented properties

The implemented assign state properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?AssignState?Default\AssignState
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?AssignState?Default\AssignState
NodeClass	Variable
BrowseName	2, "AssignState"
DisplayName	"" , "AssignState"
Description	"" , ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 18:20:39.010
SourcePicoseconds	0
ServerTimestamp	03-03-2022 18:20:39.010
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	0
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

Reference	Target Display Name
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location ($\leq \Rightarrow$ **Node/Parameter/Tag path**) in the server configuration tree:

\STATE\AssignState

4.5.2 RunState

Object values

The data type is **String**.

The defined object values (meaning the state) are as follows:

- **Ready**
- **Method Run**
- **Manual Run**
- **System Pause**
- **Hold**
- **Manual Pause**
- **Wash**
- **Alarms and errors**
- **Initializing system**
- **Resetting**
- **Starting method run**
- **Starting manual run**

The quality is good on a successful request of run state data or on a valid update from the online data, otherwise it is bad.

Implemented properties

The implemented run state properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?RunState?Default\RunState
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?RunState?Default\RunState
NodeClass	Variable
BrowseName	2, "RunState"
DisplayName	"", "RunState"
Description	"", ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	03-03-2022 18:25:39.315
SourcePicoseconds	0
ServerTimestamp	03-03-2022 18:25:39.315
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	Ready
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

\STATE\RunState

4.5.3 Command

Object values

The canonical data type is **Int32**

The defined object values are as follows:

Value	State
0	Pause
1	Hold
2	Continue
3	End
4	Next breakpoint
5	None

The method command objects are **CurrentRead** and **CurrentWrite**. Writing a valid value causes the utility object to submit a manual command.

When a client writes **END (3)**, the current run state is checked to see if the result should be saved or not. If a method is running, the run state is **PAUSE (0)** or **HOLD (1)**, and the result is saved. Other run state values imply a manual instruction is being executed, and the result is created in the [SystemName] (Manual) folder.

Implemented properties

The implemented general command properties are:

The screenshot shows the 'Attributes' window with the following data:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?Command?Default\Command
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?Command?Default\Command
NodeClass	Variable
BrowseName	2, "Command"
DisplayName	""; "Command"
Description	""; ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	04-03-2022 14:50:10.658
SourcePicoseconds	0
ServerTimestamp	04-03-2022 14:50:10.624
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	-1
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, CurrentWrite
UserAccessLevel	CurrentRead, CurrentWrite
AccessLevelEx	CurrentRead, CurrentWrite
MinimumSamplingInterval	0
Historizing	false

The 'References' section shows the following data:

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

\STATE\Command

4.5.4 InstrumentStatus

Object values

The **InstrumentStatus** handles the instrument connection state. The data type is **Int32**.

The defined object values are as follows:

Value	State
0	Ready
1	Scanning
2	Unknown

InstrumentStatus objects are **CurrentRead**.

Note: *The instrument connection must be valid (connected) for the other items to work as expected.*

Implemented properties

The implemented instrument connection properties are as follows:

Attribute	Value
NodeId	ns=2;s= 1:AvantSystem?InstrumentStatus?Default\InstrumentStatus
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?InstrumentStatus?Default\InstrumentStatus
NodeClass	Variable
BrowseName	2, "InstrumentStatus"
DisplayName	""; "InstrumentStatus"
Description	""; ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	04-03-2022 14:56:53.991
SourcePicoseconds	0
ServerTimestamp	04-03-2022 14:56:53.991
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	0
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftrightarrow **Node/Parameter/Tag path**) in the server configuration tree:

\STATE\InstrumentStatus

4.5.5 InstrumentStatusConnection

Object values

The ***InstrumentStatusConnection*** handles the instrument status state. The data type is ***Int32***.

The defined object values are as follows:

Value	Connection
0	Yes
1	Partially
2	No

InstrumentStatusConnection objects are ***CurrentRead..***

Note: *The instrument status must be valid (connected) if the other items are to work as expected.*

Implemented properties

The implemented instrument status properties are as follows:

The screenshot displays two windows from a software application. The top window, titled 'Attributes', shows a tree view of properties for a 'NodeId' and its 'Value'. The 'NodeId' properties include NamespaceIndex (2), IdentifierType (String), Identifier (1:AvantSystem?InstrumentStatusConnection?Default\InstrumentStatusConnection), NodeClass (Variable), BrowseName (2, "InstrumentStatusConnection"), DisplayName ("", "InstrumentStatusConnection"), Description ("", ""), WriteMask (0), UserWriteMask (0), RolePermissions (RolePermissionType Array[0]), UserRolePermissions (RolePermissionType Array[0]), and AccessRestrictions (None). The 'Value' properties include SourceTimestamp (04-03-2022 14:52:42.832), SourcePicoSeconds (0), ServerTimestamp (04-03-2022 14:52:42.832), ServerPicoSeconds (0), StatusCode (Good (0x00000000)), and Value (0). The 'DataType' properties include NamespaceIndex (0), IdentifierType (Numeric), Identifier (6 [Int32]), ValueRank (-1 (Scalar)), ArrayDimensions (Null), AccessLevel (CurrentRead), UserAccessLevel (CurrentRead), AccessLevelEx (CurrentRead), MinimumSamplingInterval (0), and Historizing (false). The bottom window, titled 'References', shows a table of references with columns 'Reference' and 'Target DisplayName'. The references are: HasProperty (EnumStrings) and HasTypeDefiniti... (MultiStateDiscreteType).

Attribute	Value
NodeId	ns= 2;s= 1:AvantSystem?InstrumentStatusConnection?Default\InstrumentStatusConnection
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?InstrumentStatusConnection?Default\InstrumentStatusConnection
NodeClass	Variable
BrowseName	2, "InstrumentStatusConnection"
DisplayName	"", "InstrumentStatusConnection"
Description	"", ""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	04-03-2022 14:52:42.832
SourcePicoSeconds	0
ServerTimestamp	04-03-2022 14:52:42.832
ServerPicoSeconds	0
StatusCode	Good (0x00000000)
Value	0
DataType	Int32
NamespaceIndex	0
IdentifierType	Numeric
Identifier	6 [Int32]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

\STATE\InstrumentStatusConnection

4.5.6 UserInControl

Object values

UserInControl handles the instrument control state. The canonical data type is **String**. The defined object value is the user name of the current user having control.

UserInControl objects are **CurrentRead**.

Implemented properties

The implemented instrument status properties are as follows:

Attribute	Value
NodeId	ns=2;s=1:AvantSystem?UserInControl?Default\UserInControl
NamespaceIndex	2
IdentifierType	String
Identifier	1:AvantSystem?UserInControl?Default\UserInControl
NodeClass	Variable
BrowseName	2, "UserInControl"
DisplayName	"", "UserInControl"
Description	""
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[0]
UserRolePermissions	RolePermissionType Array[0]
AccessRestrictions	None
Value	
SourceTimestamp	04-03-2022 14:58:44.927
SourcePicoseconds	0
ServerTimestamp	04-03-2022 14:58:44.927
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	Default
DataType	String
NamespaceIndex	0
IdentifierType	Numeric
Identifier	12 [String]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead
UserAccessLevel	CurrentRead
AccessLevelEx	CurrentRead
MinimumSamplingInterval	0
Historizing	false

Reference	Target DisplayName
HasProperty	EnumStrings
HasTypeDefiniti...	MultiStateDiscreteType

The following is an example location (\Leftarrow **Node/Parameter/Tag path**) in the server configuration tree:

```
\STATE\UserInControl
```

Note: **Run state** and **Assign state tag value** reflect disconnection of system once user disconnects system from OPC UA Client.

4.6 Recommendations

We recommend writing to items synchronously. The server handles both synchronous and asynchronous writes, but when making asynchronous writes, it is up to the client to wait until the server is ready with a previous asynchronous write.

- Note:**
- *The UNICORN system settings are available via the **SETTINGS** branch, which works like the **INSTRUCTION** branch with one exception. There is only one execute instruction, `\Setting\Store`, to store all system settings at once.*
 - *System settings can only be saved when the system is in **Ready** state. If UNICORN changes system settings, it updates the settings parameters in OPC.*

5 UNICORN OPC Alarms and Events address space

About this chapter

This chapter describes the UNICORN OPC Alarms and Events address space, various types of alarms, error messages, and event categories.

The Alarms and Events address space shows the internal alarm functionality of the instrument configuration. Each item in the address space might trigger an alarm or event.

Each component's failure/loss triggers an alarm and the Alarms and Events (AE) client connected in control/view of the system is notified. Alarms, events, and warnings are notified only when the user drags a **Server** object and expands or connects a node for a system under the **AlarmsEvents** folder. If the connection to the instrument is lost, then the AE client is notified through the restart error alarm. Once the restart error occurs and the connection issue has been fixed, then either the UNICORN Instrument Server computer must be restarted, or the restart error must be acknowledged in UNICORN **System Control**.

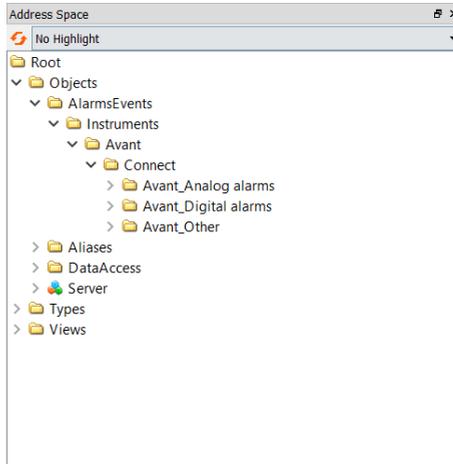
To acknowledge alarms or warnings from the OPC UA Client, the OPC UA Client (UNICORN OPC Server) must be in control of the system. Use the **Data Access** address space to take control over the system via the **AssignState** item.

The UNICORN OPC and Alarms and Events (AE) server does not support the inactive/acknowledged/enabled state. The UNICORN OPC AE server will not be able to notify that the alarm or warning becomes inactive. The UNICORN OPC AE server only notifies if the signal reaches an alarm or warning state, but it will not notify the opposite when reaching the normal state.

Note: *All alarms raised from the same source will not be available to acknowledge. Only the last alarm, warning, or error generated from the same source will be available to acknowledge. However, all the alarms and warnings will be logged under the events tab. For example, when alarms 1 to 10 are generated from the same source, only alarm 10 will be available to acknowledge. The alarms disappear from the queue once they are acknowledged.*

Alarms and Events address space view

The following screenshot shows the Alarms and Events address space view.



Tip: *The address space depends on the instrument configuration. Different items will appear depending on which instrument configuration is used.*

In this chapter

Section		See page
5.1	Alarms and errors	82
5.2	Event categories	84

5.1 Alarms and errors

Analog alarm

There are always four conditions for analog alarm objects (sources). Corresponding condition names are as follows:

1. **LO_ALARM**
2. **LO_WARNING**
3. **HI_ALARM**
4. **HI_WARNING**

Both **Condition Events** and **Tracking Events** can be generated. **Condition Events** are generated when a condition becomes active or inactive and when a condition is acknowledged. **Tracking Events** are generated when a condition is disabled but still active or inactive and unacknowledged. The event categories are **Level** for **Condition Events** and **Enable/Disable** for **Tracking Events**. No attributes are supported for analog alarms.

Example location (<=> **OPC qualified source name**) in server area space:

```
Analog alarms.Cond2
```

Digital alarm

There is only one condition for each digital alarm object (source). The condition names for the two different types are:

1. **WARNING**
2. **ALARM**

Both **Condition Events** and **Tracking Events** can be generated. **Condition Events** are generated when a condition becomes active or inactive and when a condition is acknowledged. **Tracking Events** are generated when a condition is disabled but still active or inactive and unacknowledged. Event categories are **DiscreteW** or **DiscreteA** for **Condition Events** and **Enable/Disable** for **Tracking Events**. No attributes are supported for digital alarms.

Example location (<=> **OPC qualified source name**) in server area space:

```
Digital alarms.FlowWarning
```

Errors

This is the error dialog box in UNICORN **System Control**. The event category of the condition error event is **Errors**.

No attributes are defined for error events and no acknowledgements are required for this event type.

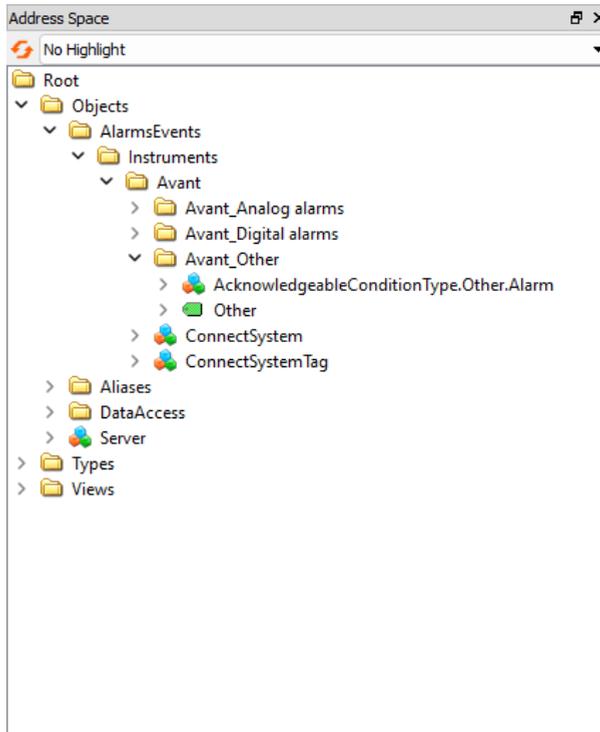
Location (<=> **OPC qualified source name**) in server area space:

Other.Errors

5.2 Event categories

Alarms and Events address space view

The following screenshot shows the Alarms and Events address space view.



6 UNICORN OPC Historical Data Access address space

About this chapter

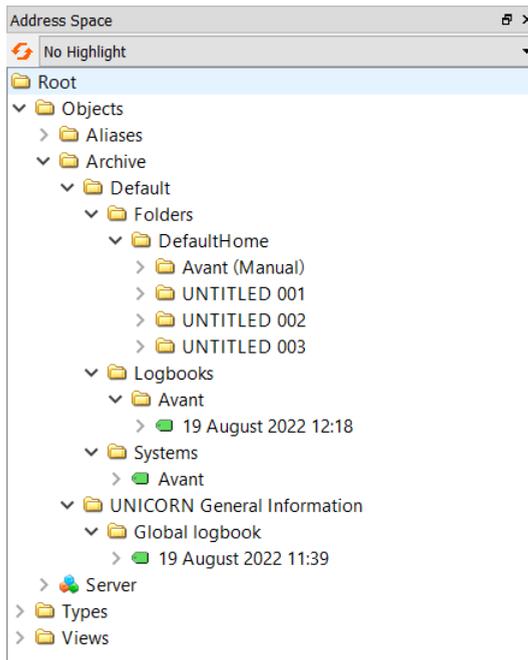
This chapter provides information about the UNICORN OPC Historical Data Access (HDA) address space, audit trail, HDA XML format definition, etc. The HDA server supports Read raw data.

The HDA address space is dynamic. A result file can be browsed whenever a new result is created. If a result is deleted, it is removed from the address space.

Tip: *The address space depends on the result files that have been created based on a specific instrument configuration. Different items will appear in the result file depending on the instrument configuration used when the result file was created.*

Historical Data Access address space view

The following screenshot shows the HDA address space view.



Note: *The address space depends on the result files that have been created based on a specific instrument configuration. Different items will appear in the result file depending on which instrument configuration that was used when the result file was created.*

In this chapter

Section		See page
6.1	AuditTrail	87
6.2	Result file information	88
6.3	UNICORN OPC Historical Data Access XML format definition	94

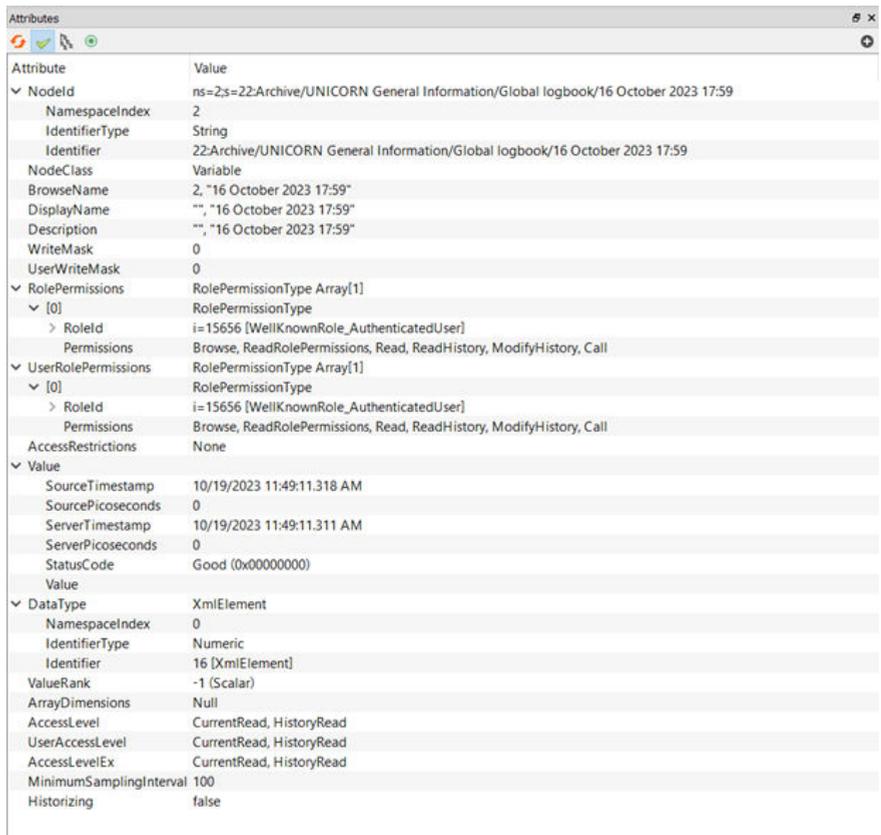
6.1 AuditTrail

Introduction

The **AuditTrail** from UNICORN is available under the HDA Server address space. The global logbook is found under the **UNICORN** → **GENERAL INFORMATION** branch and the system specific logs are found under the **Logbooks** folder in the user (e.g., **Default**) folder.

Attributes defined by **AuditTrail**

Each **AuditTrail** leaf (item) has the following attributes defined:



Attribute	Value
NodeId	ns=2s=22:Archive/UNICORN General Information/Global logbook/16 October 2023 17:59
NamespaceIndex	2
IdentifierType	String
Identifier	22:Archive/UNICORN General Information/Global logbook/16 October 2023 17:59
NodeClass	Variable
BrowseName	2, "16 October 2023 17:59"
DisplayName	"" , "16 October 2023 17:59"
Description	"" , "16 October 2023 17:59"
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[1]
[0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
UserRolePermissions	RolePermissionType Array[1]
[0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
AccessRestrictions	None
Value	
SourceTimestamp	10/19/2023 11:49:11.318 AM
SourcePicoseconds	0
ServerTimestamp	10/19/2023 11:49:11.311 AM
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
DataType	XmlElement
NamespaceIndex	0
IdentifierType	Numeric
Identifier	16 [XmlElement]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, HistoryRead
UserAccessLevel	CurrentRead, HistoryRead
AccessLevelEx	CurrentRead, HistoryRead
MinimumSamplingInterval	100
Historizing	false

The start and end time are equal for a leaf. The time is set to the creation time of the audit trail file, meaning the renew time.

6.2 Result file information

Introduction

Each result file contains **Curves**, **Documentation**, and **Peak tables**.

Information about the system is available under the **Systems** folder in the user folder.

In this section

Section	See page
6.2.1 Curves	89
6.2.2 Documentation	90
6.2.3 Peak tables	92
6.2.4 Systems	93

6.2.1 Curves

Introduction

All curves in a result file are listed as items directly under the result file branch in the address space.

Attributes supported by Curves

Each curve has several supported attributes (see below) that allow the kind of curve that the item represents, to be identified.

Attribute	Value
▼ NodeId	ns=2;s=13:Archive/Default/Folders/DefaultHome/DefaultHome/UNICORN 6.2 Avant 150/pH 3 5/pH 3 5:Chrom.1:% Cond
NamespaceIndex	2
IdentifierType	String
Identifier	13:Archive/Default/Folders/DefaultHome/DefaultHome/UNICORN 6.2 Avant 150/pH 3 5/pH 3 5:Chrom.1:% Cond
NodeClass	Variable
BrowseName	2, "pH 3 5:Chrom.1:% Cond"
DisplayName	"", "pH 3 5:Chrom.1:% Cond"
Description	"", "pH 3 5:Chrom.1:% Cond"
WriteMask	0
UserWriteMask	0
▼ RolePermissions	RolePermissionType Array[1]
▼ [0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
▼ UserRolePermissions	RolePermissionType Array[1]
▼ [0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
AccessRestrictions	None
▼ Value	
SourceTimestamp	10/19/2023 11:54:27.792 AM
SourcePicoseconds	0
ServerTimestamp	10/19/2023 11:54:27.778 AM
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
▼ DataType	Float
NamespaceIndex	0
IdentifierType	Numeric
Identifier	10 [Float]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, HistoryRead
UserAccessLevel	CurrentRead, HistoryRead
AccessLevelEx	CurrentRead, HistoryRead
MinimumSamplingInterval	100
Historizing	false

6.2.2 Documentation

Introduction

Evaluation in UNICORN has a documentation dialog for each result file. The **DOCUMENTATION** branch for each result file in the HDA contains the same data as **Evaluation** in documentation.

Items in Documentation

The following items are available:

Item	Data type
BatchNumber	String
BufferPro	XML string
Calibration	XML string
Columns	XML string
Evaluation logbook	XML string
Evaluation procedures	XML string
FracXY	XML string
Run logbook	XML string
Method name	String
Method creator	String
Method creation date	String
Method created for system	String
Method last modifier	String
Method last modification date	String
Method signatures	XML string
Method notes	XML string
Start notes	XML string
Run notes	XML string
Evaluation notes	XML string
Method strategy notes	String
Result name	String

Item	Data type
<i>Result creator</i>	<i>String</i>
<i>Result creation date</i>	<i>String</i>
<i>Result run system</i>	<i>String</i>
<i>Batch number</i>	<i>String</i>
<i>Result signatures</i>	<i>XML string</i>
<i>Result strategy used components</i>	<i>XML string</i>
<i>Scouting</i>	<i>XML string</i>
<i>System Settings</i>	<i>XML string</i>
<i>Snapshot</i>	<i>XML string</i>
<i>Text Instructions</i>	<i>XML string</i>
<i>Variables</i>	<i>XML string</i>
<i>Method Instrument Configuration</i>	<i>XML string</i>
<i>Result Instrument Configuration</i>	<i>XML string</i>
<i>Questions</i>	<i>XML string</i>

6.2.3 Peak tables

The **PEAK TABLES** branch contains all peak tables stored in the result file. When reading from the item, the peak table is returned as an XML-formatted string.

6.2.4 Systems

Introduction

Information about the system is available under the **Systems** folder in the user folder.

Items in Systems

Each system leaf defines the following attributes:

Attributes	
Attribute	Value
NodeId	ns=2;s=21:Archive/Default/Systems/Wave
NamespaceIndex	2
IdentifierType	String
Identifier	21:Archive/Default/Systems/Wave
NodeClass	Variable
BrowseName	2, "Wave"
DisplayName	"", "Wave"
Description	"", "Wave"
WriteMask	0
UserWriteMask	0
RolePermissions	RolePermissionType Array[1]
[0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
UserRolePermissions	RolePermissionType Array[1]
[0]	RolePermissionType
> RoleId	i=15656 [WellKnownRole_AuthenticatedUser]
Permissions	Browse, ReadRolePermissions, Read, ReadHistory, ModifyHistory, Call
AccessRestrictions	None
Value	
SourceTimestamp	10/19/2023 12:44:53.330 PM
SourcePicoseconds	0
ServerTimestamp	10/19/2023 12:44:53.302 PM
ServerPicoseconds	0
StatusCode	Good (0x00000000)
Value	
DataType	XmlElement
NamespaceIndex	0
IdentifierType	Numeric
Identifier	16 [XmlElement]
ValueRank	-1 (Scalar)
ArrayDimensions	Null
AccessLevel	CurrentRead, HistoryRead
UserAccessLevel	CurrentRead, HistoryRead
AccessLevelEx	CurrentRead, HistoryRead
MinimumSamplingInterval	100
Historizing	false

6.3 UNICORN OPC Historical Data Access XML format definition

Introduction

The XML format is defined for each result file item. This section describes the different XML structures.

In this section

Section	See page
6.3.1 BufferPro	96
6.3.2 Calibration	97
6.3.3 EvaluationLog	98
6.3.4 Run logbook	99
6.3.5 SignatureList	100
6.3.6 SnapshotList	101
6.3.7 UsedComponents	102
6.3.8 Notes	103
6.3.9 ScoutingList	104
6.3.10 SettingsList	105
6.3.11 TextInstructions	106
6.3.12 VariableList	107
6.3.13 QuestionList	108
6.3.14 PeakTable	109
6.3.15 Unicorn raw	112
6.3.16 Columns	114
6.3.17 EvaluationProcedures	116
6.3.18 FracXY	117
6.3.19 AuditTrail	118
6.3.20 Method/Result Instrument Configuration	119
6.3.21 System	120

6.3.1 BufferPro

Introduction

This XML structure is exported by reading the **BufferPro** leaf.

Note: For migrated UNICORN 5.x results, the **BufferPro** OPC item contains **BufferPrep** data. Both the S function and the S function share the same storage format.

BufferPro XML structure

Tag name	Description
BufferPro	Root item
RecipeName	Recipe name
pHRange	pH range, separated with "-"
B100	100% buffer B
AcidOrBase	Acid or base name
Salt	Salt name and salt stock concentration
Notes	Recipe name
Buffer	Specifications of the different buffers
Name	Buffer name
pKa1	pK_a1
pKa2	pK_a2
pKa3	pK_a3
dpKa1dt	dpk_a1 delta
dpKa2dt	dpk_a2 delta
dpKa3dt	dpk_a3 delta
AcidicProtons	Number of acidic protons
ChargeDeprotonatedIon	Number of charge deprotonated ions
SaltName	Salt name
ChargeAnion	Number of charged anions
ChargeCation	Number of charged cations

6.3.2 Calibration

Introduction

This XML structure is exported by reading the **Calibration** leaf.

Calibration XML structure

Tag name	Description
CalibrationList	Root item
Calibration	A calibration
Tagname	Name of item that has been calibrated
Date	Date
Username	User name
Point1	Point 1
Point2	Point 2
Constant	Constant
Offset	Offset
Remote	Calibration local or remote. (YES/NO)
P	Proportional control component
I	Integral control component
D	Derivative control component
ReferenceValue1	Reference value 1
ReferenceValue2	Reference value 2
CalibrationConstantDescription1	Calibration constant description 1
CalibrationConstantValue1	Calibration constant value 1
CalibrationConstantDescription2	Calibration constant description 2
CalibrationConstantValue2	Calibration constant value 2
CalibrationConstantDescription3	Calibration constant description 3
CalibrationConstantValue3	Calibration constant value 3

6.3.3 EvaluationLog

Introduction

This XML structure is exported by reading the **Evaluation logbook** leaf.

EvaluationLog XML structure

Tag name	Description
EvaluationLog	Root item.
Item	Item containing evaluation logbook event.

6.3.4 Run logbook

Introduction

This XML structure is exported by reading the **Run logbook** leaf.

Logbook XML structure

Tag name	Description
Logbook	Root item
Item	A logbook item
AccumulatedTime	Accumulated time in seconds
AccumulatedVolume	Accumulated volume, including unit
Event	Logbook event
EventType	Event type
EventSubType	Subtype of event

6.3.5 SignatureList

Introduction

This XML structure is exported by reading the **Method Signatures** and **Result Signatures** leaves.

SignatureList XML structure

Tag name	Description
SignatureList	Root item
Signature	A signature
Username	Name of user
Fullname	Full name of user
Position	Position of user
Date	Date, formatted in server local time
Meaning	Signature meaning
Locked	If present, the file is locked against further change

6.3.6 SnapshotList

Introduction

This XML structure is exported by reading the **Snapshot** leaf.

Note: *To get the actual y-axis value, use the time or volume value as a reference into the curves.*

SnapshotList XML structure

Tag name	Description
SnapshotList	Root item
Chromatogram	A signature
Name	Name of chromatogram
Snapshot	Run number
TimeRetention	Time retention
VolumeRetention	Volume retention
Curve	Curve
Name	Curve name
Unit	Curve unit
Time Value	Time stamp of snapshot
Volume Value	Volume stamp of snapshot

6.3.7 UsedComponents

Introduction

This XML structure is exported by reading the **Result Strategy Used Components** leaf.

UsedComponents XML structure

Tag name	Description
<i>UsedComponents</i>	Root item.
<i>Component</i>	Component name used during run.

6.3.8 Notes

Introduction

This XML structure is exported by reading the **Method Notes**, **Start Notes**, **Run Notes**, and **Evaluation Notes** leaves.

Notes XML structure

Tag name	Description
Notes	Root item, the note follows.

6.3.9 ScoutingList

Introduction

This XML structure is exported by reading the **Scouting** leaf.

Note: *To get the actual y-axis value, use the time or volume value as a reference into the curves.*

ScoutingList XML structure

Tag name	Description
ScoutingList	Root item.
Scouting	A scouting run.
TotalNumberOfScoutings	Total number of scouting items available.
RunScouting	If available, this item indicates the scouting run number of this result file.
Run	Run number.
ThisScoutingWasUsedDuringRun	This leaf is available if this scouting run was used when the method was run. The run number is the same as RunScouting .
Variable	Variables defined in this run.
Block	Block name of this variable.
Name	Name of variable.
Unit	Unit, only included if variable has a unit.
Value	Value of variable.

6.3.10 SettingsList

Introduction

This XML structure is exported by reading the **System Settings** leaf.

SettingsList XML structure

Tag name	Description
SettingsList	Root item.
Group	A group.
GroupName	The name of the group.
Instruction	Instruction names. There is at least one instruction per group.

6.3.11 TextInstructions

Introduction

This XML structure is exported by reading the ***Text Instructions*** leaf.

TextInstructions XML structure

Tag name	Description
<i>TextMethod</i>	Root item, the complete text method follows. Each row is separated by a carriage return and a new line.

6.3.12 VariableList

Introduction

This XML structure is exported by reading the **Variables** leaf.

Start protocol variables XML structure

Tag name	Description
VariableList	Root item.
Variable	A variable.
Block	Block name of the variable.
Name	The variable name.
VisibleInScouting	Indicates if the variable is visible in scouting.
VisibleInDetails	Indicates if the variable is visible in details only.
Value	Value. Can be a number or a string, depending on the variable.
Unit	Unit of the value. Not included if no unit is present.

6.3.13 QuestionList

Introduction

This XML structure is exported by reading the **Questions** leaf.

Start protocol questions XML structure

Tag name	Description
QuestionList	Root item
Item	A question item
Question	The question
Answer	The answer

6.3.14 PeakTable

Introduction

This XML structure is exported by reading leaves under the **PEAK TABLES** branch.

PeakTable XML structure

Tag name	Description
PeakTable	Root item
Summary	Summary of the peak table
TotalNumberOfDetected-Peaks	Total number of peaks
TotalArea	Total area
Unit	Unit used.
AreaInEvaluatedPeaks	Area in evaluated peaks
RationPeakareaTotalarea	Ratio peak area/total area
TotalPeakDuration	Total peak duration
ColumnHeight	Column height
ColumnV0	Column V_0
ColumnVt	Column V_t
SourceCurve	Source curve name
Baseline	Baseline
Rejection	Rejection
MinHeight	Minimum height
MinWidth	Minimum width
MaxWidth	Maximum width
MinArea	Minimum area
MaxNumberOfPeaks	Maximum number of peaks
Quantitation	Quantitation (molecular size, standard addition, internal standard, external standard, recovery)
Peak	A peak

Tag name	Description
No	Peak number
Name	Peak name
Retention	Retention
Start	Start
End	End
Width	Width
Area	Area
AreaPerTotalArea	Area/total area
AreaPerPeakArea	Area/peak area
Height	Height
WidthAtHalfHeight	Width at half peak height
HeightAtStart	Height at start
HeightAtEnd	Height at end
BaselineAtStart	Baseline at start
BaselineAtMax	Baseline at max
BaselineAtEnd	Baseline at end
ConductivityHeightStart	Conductivity height at start
ConductivityHeightMax	Conductivity height at maximum
ConductivityHeightEnd	Conductivity height at end
FractionTubeAtStart	Fraction tube at start
FractionTubeAtMax	Fraction tube at maximum
FractionTubeAtEnd	Fraction tube at end
PeakStartType	Peak start type (dropline, skim, baseline, skim-drop, unknown)
PeakEndType	Peak end type (dropline, skim, baseline, skim-drop, unknown)
Sigma	Sigma
Resolution	Resolution
CapacityFactor	Capacity factor

Tag name	Description
<i>Kav</i>	K_{av}
<i>PlateHeight</i>	Plate height
<i>PlatesPerMeter</i>	Plates per meter
<i>AsymmetryStart</i>	Asymmetry start
<i>AsymmetryEnd</i>	Asymmetry end
<i>Asymmetry</i>	Asymmetry
<i>AverageConductivity</i>	Average conductivity
<i>ExtinctionCoefficient</i>	Extinction coefficient
<i>ExtCoeffConcentration</i>	Concentration calculated using extinction coefficient.
<i>ExtCoeffAmount</i>	Amount calculated using extinction coefficient.
<i>Concentration</i>	Concentration
<i>Amount</i>	Amount
<i>MolecularSize</i>	Molecular size
<i>Type</i>	Type

6.3.15 Unicorn raw

Introduction

This XML structure is exported by reading the attribute **HDA_UNICORN_RAW_DATA** for a **Curve** leaf.

UnicornRawData XML structure

Tag name	Description
UnicornRawData	Root item.
Name	Name of curve.
Type	Curve type. Does not contain the bit mask, which the HDA_CURVE_TYPE does, only the actual type.
Unit	Curve unit (y-axis).
TimeUnit	Unit of time curve (x-axis).
VolumeUnit	Unit of volume curve (x-axis).
CurveMin	Minimum value of y-axis.
CurveMax	Maximum value of y-axis.
ZeroAdjust	Present if zero adjust is enabled.
ZeroAdjustTime	Zero adjust time.
ZeroAdjustVolume	Zero adjust volume.
StartTimeRetention	Retention start time.
InjectionTimeRetention	Injection start time.
NumberOfTimeDataPoints	Number of curve time data points.
CTD	Curve Time Data. This is repeated for all time data points. The values are assigned as attributes to the tag (x and y attributes).
StartVolumeRetention	Retention start volume.
InjectionVolumeRetention	Injection start volume.

Tag name	Description
NumberOfVolumeDataPoints	Number of curve volume data points.
ColumnVolume-Defined	Present if column volume is defined.
ColumnVolume	Column volume value for curve.
ColumnVolumeUnit	Unit of ColumnVolume curve (X-axis).
CVD	Curve Volume Data. This is repeated for all volume data points. The values are assigned as attributes to the tag (x , y and, if available, cv).
CMD	Curve Markers Data. This is repeated for all marker (text) data points. The values are assigned as attributes to the tag (t = time, v = volume, t1 = text1 and, if available, t2 = text2).

6.3.16 Columns

Introduction

This XML structure is exported by reading the **Columns** leaf.

Columns XML structure

Tag name	Description
Columns	Root item.
Column	A column.
Name	Name of column.
Height	Height of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
Diameter	Diameter of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
Volume	Volume of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
VolumeUnit	Height of column. Attribute mandatory defines if item is mandatory.
Technique	Technique used by column. Attribute mandatory defines if item is mandatory.
Vt	Total volume of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
Vo	Empty volume of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
MaxPressure	Maximum pressure of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
DefaultFlowrate	Default flow rate of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
MaxFlowrate	Maximum flow rate of column. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
TypPeakWidthAtBase	Typical peak width at base. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
pHLongMax	Long-term maximum pH. Attribute mandatory defines if item is mandatory.

Tag name	Description
<i>pHLongMin</i>	Long-term minimum pH. Attribute mandatory defines if item is mandatory.
<i>pHShortMax</i>	Short-term maximum pH. Attribute mandatory defines if item is mandatory.
<i>pHShortMin</i>	Short-term minimum pH. Attribute mandatory defines if item is mandatory.
<i>AverageParticle-Diameter</i>	Average particle diameter. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.
<i>Code</i>	Code of column. Attribute mandatory defines if item is mandatory.
<i>TypicalLoading-Range</i>	Typical loading range. Attribute mandatory defines if item is mandatory.
<i>MolWeight-Range</i>	Molecular weight range. Attribute mandatory defines if item is mandatory. Attribute unit defines unit of value.

6.3.17 EvaluationProcedures

Introduction

This XML structure is exported by reading the **Evaluation procedure** leaf.

EvaluationProcedures XML structure

Tagname	Description
EvaluationProcedures	Root item.
EvaluationProcedure	An evaluation procedure.
Name	Name of evaluation procedure.
WasRun	Indicates if the evaluation procedure was run during method run.

6.3.18 FracXY

Introduction

This XML structure is exported by reading the **FracXY** leaf.

*FracXY*XML structure

Tagname	Description
<i>FracXY</i>	Root item
<i>FractionOrder</i>	Fraction order (Serpentine row, Row-by-row, Serpentine column or Column-bycolumn)
<i>Name</i>	Name of rack
<i>Group</i>	Rack group
<i>Name</i>	Name of rack group
<i>LastTube</i>	Last tube of group

6.3.19 AuditTrail

Introduction

This XML structure is exported by reading a **Global Logbook** or **System Logbook** leaf.

AuditTrail XML structure

Tag name	Description
AuditTrail	Root item.
Type	Type of audit trail: Global , System or Backup . Backup is used when accessing a direct audit trail file address.
ChecksumError	If this item is available, the file has been modified outside UNICORN or is corrupt.
Audit	An audit item.
Time	Time of audit.
TimeZone	Time zone.
Message	Audit message.
Details	Log entry details.

6.3.20 Method/Result Instrument Configuration

Introduction

This XML structure is exported by reading a **Method Instrument Configuration** or **Result Instrument Configuration** leaf.

Method/Result Instrument Configuration XML structure

Tag name	Description
InstrumentConfiguration	Root item
Name	Name of instrument configuration
Version	Instrument configuration version
Help text	Instrument configuration help text
StrategyName	The instrument configuration name
StrategyNotes	The instrument configuration notes
PhaseConfiguration	Phase configuration name and version
TemplateMethodConfiguration	Protocol configuration name and version
PerformanceTestConfiguration	Performance test configuration name and version
FlowSchemeConfiguration	Flow scheme configuration name and version
VIDConfiguration	VID configuration name and version
FirmwareConfiguration	Firmware configuration name and version
CompatibleUNICORNClientVersion	UNICORN version
CompatibleUNICORNInstrumentServerVersion	UNICORN version

6.3.21 System

Introduction

This XML structure is exported by reading a **System** leaf.

System XML structure

Tag name	Description
SystemName	Name of system.
SystemType	Type of system.
InstrumentConfiguration	Instrument configuration name and version.
SystemIsNonActive/SystemIsActive	Active system.
ControlledByUNICORNInstrumentServer	Instrument server computer name. Note: <i>Displayed only if the system is Active.</i>
InstrumentSerialNo	Instrument serial number.
ConnectionBasedOnSerial/ConnectionBasedOnIP	Type of connection based on serial number or instrument IP address.
InstrumentIPAddress	Instrument IP address.

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 Corporation or an affiliate doing business as Cytiva.

ÄKTA, ÄKTA process, ReadyToProcess, and UNICORN are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Microsoft and Windows are trademarks of the Microsoft group of companies.

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

© 2024 Cytiva

UNICORN© 2024 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

29752660 AA V:1 05/2024