



Automation Design Document for the Agilent Seahorse XF Pro Analyzer



Notices

Publication Number

5994-4689EN Rev A

February 25, 2022

Copyright

© Agilent Technologies, Inc. 2022

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

Contact Information

Agilent Technologies, Inc.
5301 Stevens Creek Boulevard
Santa Clara, CA 95051
USA

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

Contact page:
<https://www.agilent.com/en/contact-us/page>

Documentation feedback:
documentation.automation@agilent.com

Acknowledgements

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

Warranty

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

Technology Licenses

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

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data -Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

WARNING

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

CAUTION

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

Contents

- 1. Purpose 1
- 2. Using the Automation Client to connect to the XF Pro Controller 3
- 3. Automation Service and XF Pro Controller software 7
- Automation Properties 8
- Automation Methods 9
- 4. Error Codes 13

This page intentionally blank

1 Purpose

The purpose of this document is to define an interface between the XF Pro Controller software on a host XF Pro Controller (the computer connected to the XF Pro Analyzer) and your custom-built or third-party automation client software, for purposes of automating your XF data acquisition workflow. The host service will run on XF Pro Controller software on an XF Pro Analyzer, and the client is designed to be run separately.

The scheduling software is responsible for calling the server methods and coordinating robot actions for XF Pro cell plate and sensor cartridge handling. A valid XF Discovery software license is required to unlock the automation service for the XF Pro Analyzer. This automation service is designed for operation with one client at a time.

NOTE

Although this document references the term “XFE” throughout, the automation service is supported on XF Pro Analyzers and XF Pro Controllers only. No other Agilent Seahorse Analyzers feature an automation service as described in this document.

1 Purpose

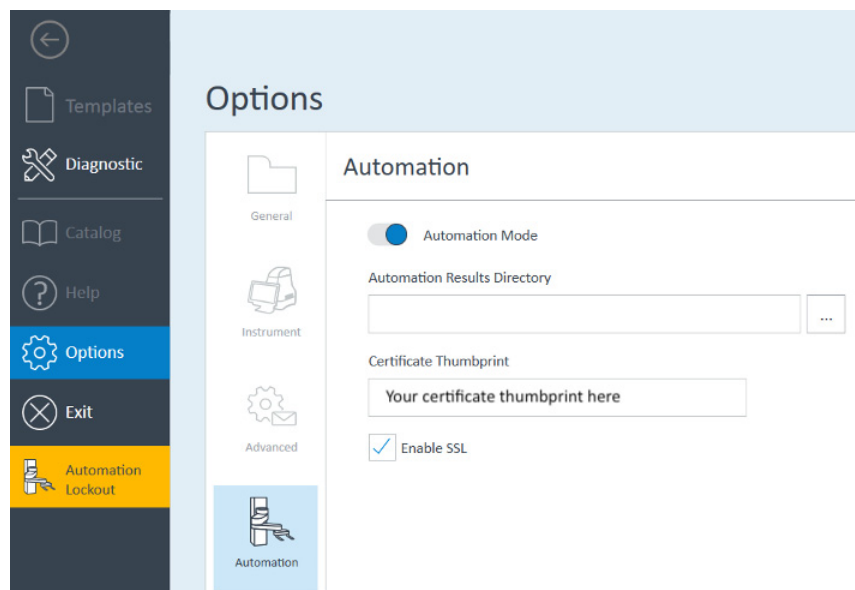
This page intentionally blank

2 Using the Automation Client to connect to the XF Pro Controller

Only an XF Pro Analyzer + Controller with a valid XF Discovery license can access the automation service. The XF Pro Controller software must be run in administration mode to run the automation service.

To enable automation mode:

- 1 Open the XF Pro Controller software.
- 2 Select **Options > Automation**.
- 3 Enable **Automation Mode**.
- 4 Enter the **Certificate Thumbprint** field and **Enable SSL**.
- 5 To apply the changes, select **Save**. Onscreen buttons are now disabled and **Automation Lockout** is displayed.



After configuring the automation service in XF Pro Controller software, you may encounter a Windows Firewall warning when starting up the Automation service (see [Figure 1](#)). Ensure the check boxes for *Domain networks* and *Private networks* are selected and click **Allow access**. This will allow XF Pro Controller software to access port 8001 and Windows Firewall will automatically create a new rule for it to access the port.

2 Using the Automation Client to connect to the XF Pro Controller

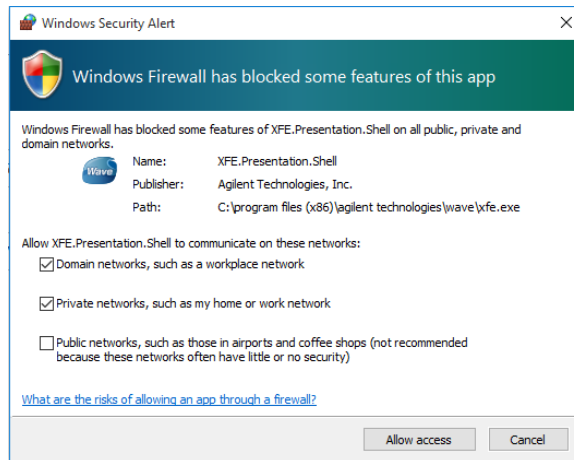


Figure 1. Windows Firewall warning.

To verify this worked successfully, open a web browser (IE recommended) and go to: [https://\[CN\]:8001/CellAnalysisServices/](https://[CN]:8001/CellAnalysisServices/)

Example: [https://\[server computer name\].\[sub domain\].\[domain\]:8001/CellAnalysisServices/](https://[server computer name].[sub domain].[domain]:8001/CellAnalysisServices/) where from the certificate, CN=[server computer name].[sub domain].[domain]

If successful, it should show up in the URL as shown in **Figure 2**.

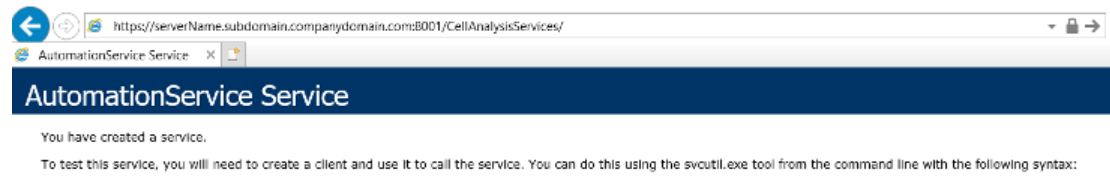


Figure 2. Successful Service.

NOTE

To test the connection, Internet Explorer is recommended. Chrome or Edge may display an erroneous warning that the connection is not private because the certificate uses a common name rather than a Subject Alternative Name (SAN) (see Figure 3).

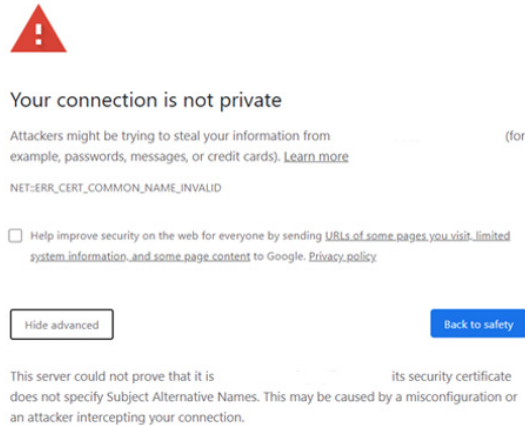


Figure 3. Browser connection warning.

Once automation mode is enabled, XF Pro Controller software is ready to begin receiving method calls from a client application interface. The client application running on a server or separate laptop or computer communicates with the XF Pro Controller through WsHTTP protocol.

On a local network, the IP address and port is needed of the XF Pro Controller in order for the client application to connect. To create an instance of the XFeAutomationClient class with the IP address and port as parameters:

```
AutomationClient client = new AutomationClient("111.111.111.111", 8000);
```

In the C# example, the IP address of the host controller is 111.111.111.111 and the port is 8000. If you are using SSL, use port 8001. The IP Address should be input as a string, following standard IP address formatting, and the port as an integer.

2 Using the Automation Client to connect to the XF Pro Controller

This page intentionally blank

3 Automation Service and XF Pro Controller software

The automation service serves as an interface to XF Pro Controller software. The interface to the automation service is composed of several methods that allow for automating data acquisition on the XF Pro Analyzer. The automation client application allows access to these methods which in turn control the XF Pro Analyzer and Controller software while the automation service is running. There are several properties of the automation service that represent basic statuses of the XF Pro Analyzer + Controller.

Automation Properties

XFeState

Parameter	Value
Type	String
Access	Read Only
Description	The current state of the XF Pro Analyzer
Example	"Ready"

Tray Status

Parameter	Value
Type	String
Access	Read Only
Description	The current state of the XF Pro Tray
Example	"In"

Automation Methods

GetStatus

The GetStatus method returns the current state of the instrument.

```
string GetStatus()
```

Return value: The instrument status

XFeState Enumeration	String Equivalent	Meaning
Ready	Ready	The instrument is ready to run an assay.
Running	Running	The instrument is running an assay.
RunningWaitingForReadyToLoadCartridgeAndCalibrantPlate	Ready to Load Cartridge and Plate	The instrument is ready for the user to open the tray for the cartridge and calibration plate to be loaded.
RunningWaitingForLoadCartridgeAndSupportingPlate	Waiting for Cartridge and Plate	The tray is open, and the instrument is ready for the user to place the cartridge and calibration plate on the tray.
RunningLoadCartridgeAndCalibrantPlate	Loading Cartridge and Plate	The user has loaded the cartridge and calibrant plate on the tray and has called LoadCartridge.
RunningCalibration	Calibrating	The instrument is running Calibration.
RunningCalibrationComplete	Finished Calibration	The instrument has finished Calibration.
RunningWaitingForReadyToLoadCellPlate	Ready to Load Cell Plate	The instrument is ready for the user to open the tray for the cell plate to be loaded.
RunningWaitingForLoadCellPlate	Waiting for Cell Plate	The tray is open, and the instrument is ready for the user to place cell plate on the tray.
RunningLoadCellPlate	Running	The instrument is running the Assay.
RunningComplete	Assay Finished	The Assay has completed successfully.
RunningWaitingForReadyToStripCartridge	Ready to Unload Cartridge	Assay has finished (or cancelled), ready to strip/unload cartridge message pops up.
RunningWaitingForStripCartridge	Waiting for Unload Cartridge	The instrument is waiting for the user to unload the cartridge.
RunningStripCartridge	Unloading Cartridge and Plate	Cartridge and Cell plate have been stripped and unloaded from tray, assay has completed.
ReadyRunComplete	Assay Completed!	The Assay has finished and is about to transition back to the Ready state.
AssayCancelled	Assay Cancelled!	The user has canceled the assay. The instrument will transition back to Ready state.

GetTrayStatus

The GetTrayStatus method returns the current state of the tray.

string GetTrayStatus()

Return value: The tray status

TrayStatus Enumeration	String Equivelant	Meaning
In	In	The tray is in.
Out	Out	The tray is out.

GetErrors

The GetErrors method returns the current list of errors.

string GetErrors()

Return value: Error text

OpenTray

The OpenTray method opens the tray in order to load/unload the cartridge or to load the cell plate.

bool OpenTray()

Return value: Boolean if client OpenTray method call succeeds

LoadCartridge

The LoadCartridge method loads the cartridge and utility plate after an assay has started and the tray has been opened.

bool LoadCartridge()

Return value: Boolean if client LoadCartridge method call succeeds

UnloadCartridge

The UnloadCartridge method unloads the cartridge and utility plate after an assay has completed and the tray has been opened.

bool UnloadCartridge()

Return value: Boolean if client UnloadCartridge method call succeeds

LoadCellPlate

The LoadCellPlate method loads the cell plate after calibration has finished and the utility plate has been removed with the tray open.

bool LoadCellPlate()

Return value: Boolean if client LoadCellPlate method call succeeds

LoadTemplate

The LoadTemplate method loads the given XF Pro template given the input name of the template file that is saved on the controller.

bool LoadTemplate(**string** name)

Return value: Boolean if client LoadTemplate method call succeeds

Run

The Run method starts and runs the assay given the loaded template.

void Run()

Return value: Void

This page intentionally blank

4 Error Codes

The following error codes can be displayed in the error text message returned by the GetErrors client method. Most errors result in the stoppage of Automation mode and will present user-facing dialogs in XF Pro Controller software that a user must physically address to resolve the issue. If you encounter error codes not described in this list, please contact Agilent Technologies Cell Analysis Support.

Code	Label	Meaning
1	<code>XFEStripCartridgeResultType.NoCommunicationError</code>	Communication error with instrument while stripping the cartridge.
2	<code>XFEStripCartridgeResultType.GripperReleaseError</code>	The grippers did not release while stripping the cartridge.
3	<code>XFEStripCartridgeResultType.StripAborted</code>	Strip cartridge failed going to the Home position.
4	<code>XFEStripCartridgeResultType.UnknownError</code>	Unknown error while attempting to strip the cartridge.
5	<code>ControllerBase.ExceptionCaughtInMProbes</code>	Error while moving probes.
7	<code>XFEApplicationExceptionType.InjectFailure</code>	Inject failure.
8	<code>ControllerBase.ExecuteSetAgitateStatus; Value of endoperation</code>	General error during mix operation.
10	<code>ControllerBase.Calibrate</code>	Error or abort called during calibration when moving tray to the measure position.
11	<code>ControllerBase.Calibrate</code>	Error or abort called during calibration when moving probes to the Z measure position.
12	<code>ControllerBase.Calibrate</code>	Autozero failure.
13	<code>ControllerBase.Calibrate</code>	Error while attempting to run Autozero.
14	<code>ControllerBase.Calibrate; AutoLED returned an error</code>	Coarse calibration failure.
15	<code>ControllerBase.Calibrate; AutoLED Error thrown</code>	Error while attempting to run Coarse Calibration.
16	<code>XFeControllerBase.Calibrate</code>	Error while attempting to run Fine Calibration or Measure.
27	<code>XFELoadCartridgeResultType.MovementError</code>	Load cartridge failure or load plate failure or moving probes operation was aborted.
40	<code>XFELoadCartridgeResultType.LoadAborted</code>	Load cartridge failed going to the Home position.
41	<code>ControllerXFe_96.MeasureMultiBoard</code>	Measure failure.
42	<code>XFeControllerBase.ReadCartridgeBarcode</code>	No barcode reader found when loading cartridge.

Code	Label	Meaning
43	XFeControllerBase.ReadCartridgeBarcode	No cartridge barcode error when loading cartridge, or cartridge type is not available or cartridge is not compatible with instrument.
44	XFeControllerBase.ReadCartridgeBarcode	Cartridge is not compatible with the Injections in the assay protocol.
45	ControllerXFe_96.LoadCartridge	Load cartridge failed going to the Measure position.
46	XFELoadCartridgeResultType.GripperReleaseError	The grippers did not release while loading the cartridge.
47	ControllerXFe_96.LoadCartridge	Load cartridge failure while going to the load/unload position (probe head).
48	XFELoadCartridgeResultType.GripperEngageError	The grippers did not engage while loading the cartridge.
49	XFELoadCartridgeResultType.MovementError	Error while moving to the Home Z position during load cartridge.
50	XFELoadCartridgeResultType.NoCartridgeDetectedError	No cartridge detected.
51	XFELoadCartridgeResultType.LoadAborted	Load cartridge failure while going to the load position.
52	XFELoadCartridgeResultType.MonitoringError	Unable to monitor tray movement while loading cartridge.
53	ControllerXFe_96.LoadCartridge	Load Cartridge Failure, or load failure while attempting move tray to barcode position to read the plate barcode or load aborted.
54	XFELoadPlateResultType.NoBarcodeReaderError	No barcode reader detected while attempting to read the plate barcode.
55	XFELoadPlateResultType.NoBarcodeError	No plate barcode detected while attempting to read the plate barcode.
56	XFELoadPlateResultType.PlateCompatibilityError	The plate is not compatible with the instrument.
57	XFELoadPlateResultType.MonitoringError	Unable to monitor tray movement while loading plate.
58	ControllerBase.LoadPlate	Communication error while loading plate.
60	XFESTripCartridgeResultType.StripAborted	Error moving tray to measure position while attempting to strip the cartridge.
61	XFESTripCartridgeResultType.StripAborted	Error moving probes to Z home position while attempting to strip the cartridge.
64	XFESTripCartridgeResultType.StripAborted	The strip procedure was aborted while attempting to strip the cartridge.
66	ControllerBase.Measure	Error or abort called during measure when moving tray to the measure position.
67	ControllerBase.Measure	Error or abort called during measure when moving probes to the Z measure position.
68	ControllerBase.Inject	Error while moving tray to measure position during an injection.

Code	Label	Meaning
69	<code>ControllerBase.Inject</code>	Error while moving probes to inject position during an injection.
70	<code>XFEApplicationExceptionType.MixFailure</code>	Mix was aborted or error while attempting to mix.
71	<code>XFEApplicationExceptionType.MixFailure</code>	Mix error or mix was aborted while attempting to move probes to Z mix position.
80	<code>ControllerBase.Wait</code>	Error or abort called during measure wait cycle when moving tray to the measure position.
81	<code>ControllerBase.Wait</code>	Error or abort called during measure wait cycle when moving probes to the Z wait position.
83	<code>XFELoadPlateResultType.CartridgeAlreadyLoaded</code>	Cartridge detected while attempting to load plate.
97	<code>Load Plate Failure</code>	Load plate error while doing load cartridge operation.
98	<code>XFELoadCartridgeResultType.NoBarcodeError</code>	Cartridge barcode read failure.
99	<code>XFEStripCartridgeResultType.CartridgeDetectedError</code>	A cartridge was still detected after attempting to strip the cartridge.
100	<code>XFEStripCartridgeResultType.StripOK</code>	Unload cartridge failure.
101	<code>ExecuteEjectPlate</code>	Abort called while ejecting plate during unload cartridge operation.
102	<code>ExecuteEjectPlate</code>	Abort called while moving tray to home position during eject plate operation.
197	<code>XFELoadCartridgeResultType.MovementError</code>	Load Plate Failure while loading the cartridge.
311	<code>XFELoadCartridgeResultType.CartridgeAlreadyLoaded</code>	Cartridge was already loaded.

This page intentionally blank

In This Guide

This guide provides automation design interface documentation for the Agilent Seahorse XF Pro Analyzer and XF Pro Controller software including:

- Using the Automation Client to connect to the XF Pro Controller
- Automation Service and XF Pro Controller software
- Error Codes

www.agilent.com

© Agilent Technologies, Inc. 2022

February 25, 2022



5994-4689EN Revision A

