



Test Services

Administration Guide

Notices

Document Identification

Document No: D0015137
April 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.

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

Software Revision

This guide is valid for the 3.5 revision or higher of the Test Services program until superseded.

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

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.

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.

In This Guide

This guide provides administrative procedures to ensure that the Agilent Test Services tool remains operational and performs well over time. It is intended for the system administrator of the Test Services tool. Basic administrative knowledge of the OpenLab CDS software is required.

1 Introduction

This chapter describes Test Services and the QualA licensing that enables QualA tests.

2 Administration

This chapter contains information on administrative tasks for Test Services.

3 Troubleshooting

This chapter contains information on how to troubleshoot issues and resolve error messages that occur while using Test Services.

Content

1	Introduction	6
	Overview	7
	Test Services System Architecture	8
	Starting Test Services	11
2	Administration	12
	Installed Services	13
	Licensing	14
	Setting User Account to Run Test Services	15
	Test Services user roles and privileges	15
	Required roles and privileges	16
	Test Services user role creation in OpenLab ECM 3.x	17
	Logs	19
	Viewing the Test Services logs	19
	Configuring Test Services	20
	Modifying the port number	20
	If OpenLab CDS configuration changes after installation of Test Services	22
	If authentication changes after installation of Test Services	22
	Location of Test Services Test Results	23
	OpenLab Server and ECM XT	23
	ECM 3.5 and 3.6	23
	Security and Certificates	24
	Configure Test Services to use a third party certificate	24
3	Troubleshooting	26
	Troubleshooting	27
	Results reported as not passed	27

Sequence hangs during workflow test	28
Test Services projects and instruments	29
Email notification not working	29
Error Messages	31
Test Services error messages	31
OpenLab CDS Workflow Test error messages	34



1

Introduction

Overview 7

Test Services System Architecture 8

This chapter describes Test Services and the QualA licensing that enables QualA tests.

Test Services is installed automatically when you install OpenLab CDS. It includes a suite of verification tests that are available without a license. QualA for OpenLab CDS and QualA for OpenLab Server licenses enable additional CDS verification tests for the Test Services version currently installed.

For information on how to use Test Services, see the *Test Services User Guide*.

Overview

Test Services is installed when you install OpenLab CDS. It provides a framework for installation and running verification tests. Some tests are installed and enabled by default without a license. The installed Test Services CDS Plug-in contains tests that are specific to OpenLab CDS. Some of these tests require a QualA license in order to run them.

For information on how to install licenses, see [“Licensing”](#) on page 14.

NOTE

All applicable licenses are included with OpenLab CDS Workstation Plus systems. For other CDS products, licenses must be purchased and installed in order to run licensed tests.

Test Services System Architecture

Examples of Test Services deployment on CDS Workstation and Workstation Plus are shown in **Figure 1**. Examples of Test Services deployment in an OpenLab CDS Client/Server environment with OpenLab Server/ECM XT are shown in **Figure 2** and **Figure 3**. An example of Test Services for ECM storage is shown in **Figure 4**.

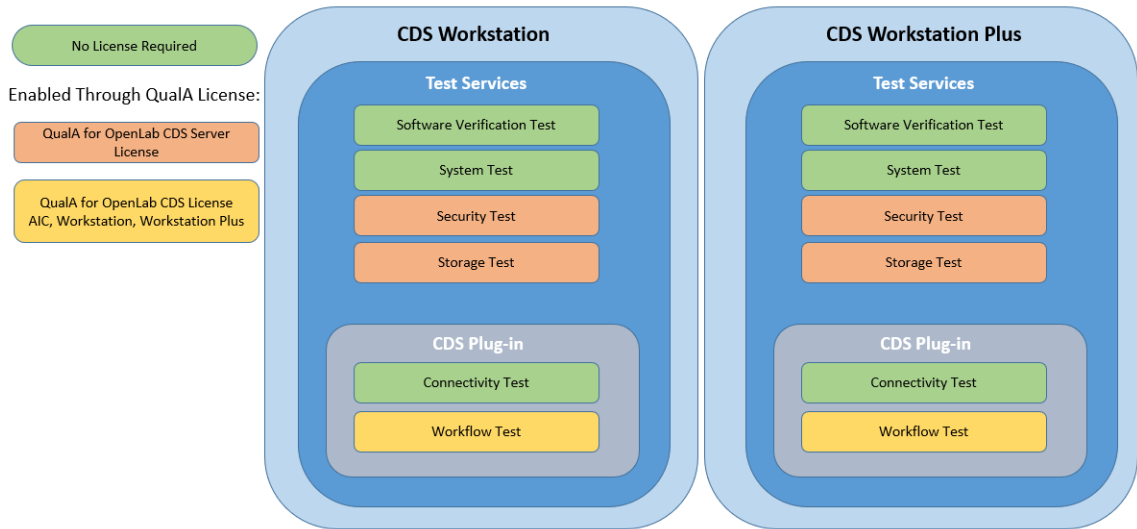


Figure 1. Test Services on Workstation or Workstation Plus

NOTE

OpenLab CDS Workstation Plus includes all QualIA licenses. For all other systems, you must purchase QualIA licenses separately.

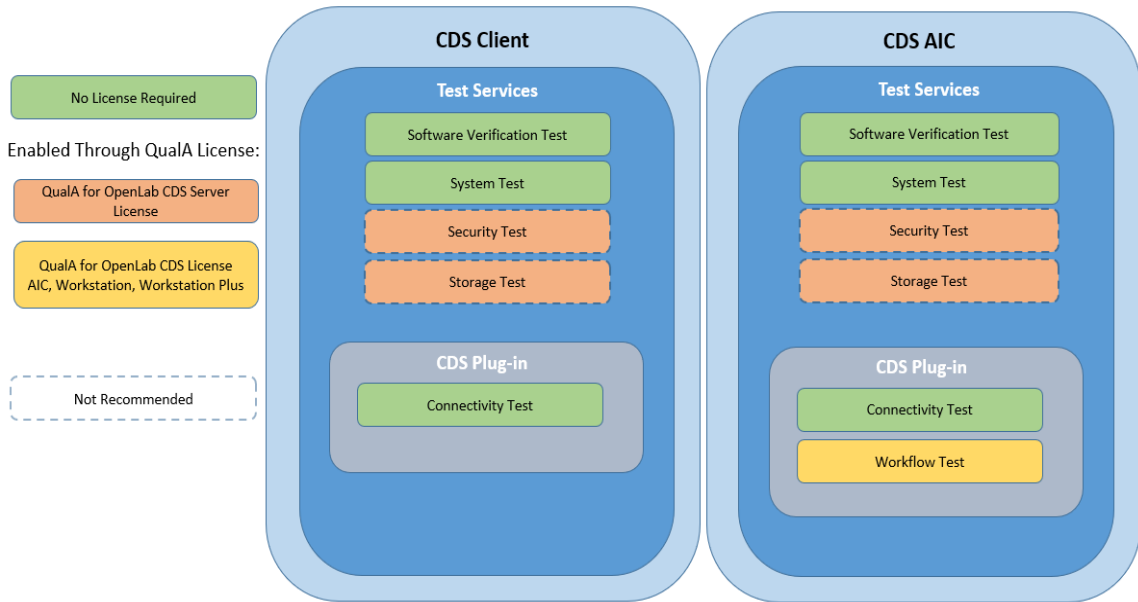


Figure 2. Test Services on CDS Client and AIC

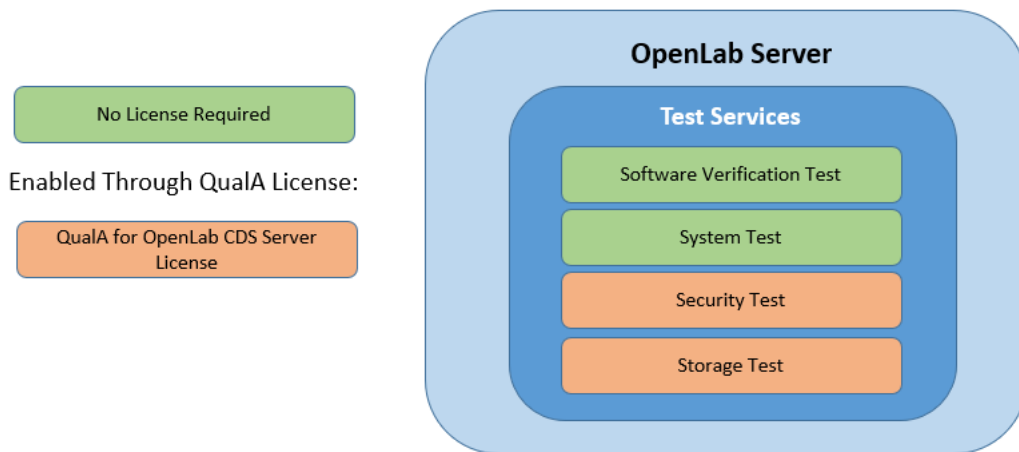


Figure 3. Test Services on OpenLab Server

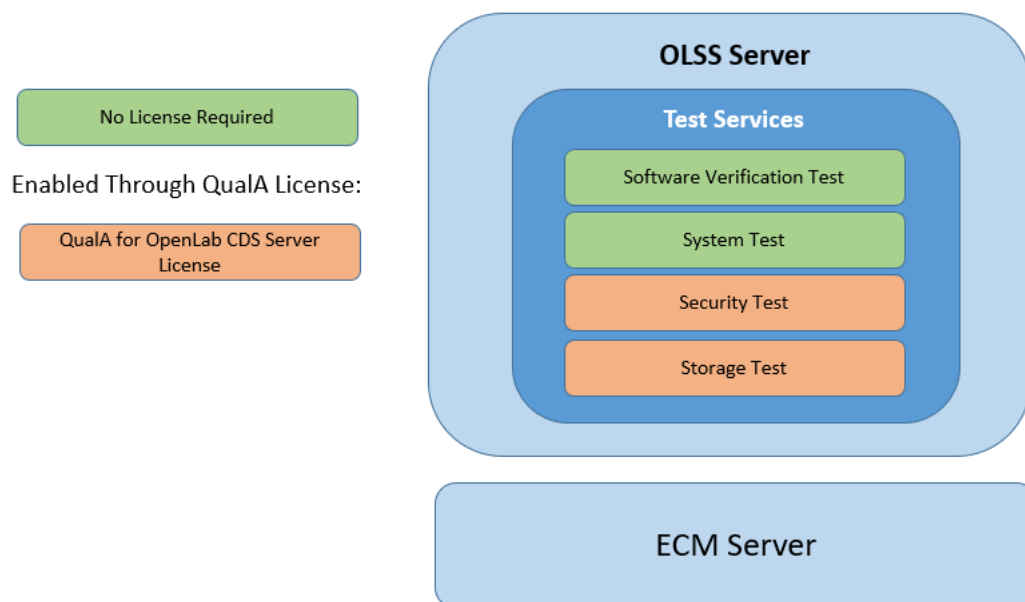


Figure 4. Test Services on OpenLab ECM

Starting Test Services

Test Services is accessed using a browser, by visiting a URL containing the machine on which Test Services is installed.

An example of the URL for opening the user interface of Test Services is shown below:

`https://<<hostname>>/testservices/`

For example: `https://localhost/testservices/`

On client systems, you must use `https://localhost:9092/` to access test services locally. If you need to access test services remotely, use `https://<<client-fqdn>>:9092/`. If there is a firewall on the client system, make sure that port 9092 is open.



2

Administration

Installed Services	13
Licensing	14
Setting User Account to Run Test Services	15
Logs	19
Configuring Test Services	20
Location of Test Services Test Results	23
Security and Certificates	24

This chapter contains information on administrative tasks for Test Services.

Installed Services

The Test Services installation creates a Windows Service named Agilent Test Services.

For OpenLab Workstation, Workstation Plus, and Server, Agilent Test Services Central Management service is also created. Agilent Test Services service installed on the Server depends on this service.

To verify the operation of the service, open Windows Run, type `services.msc`, and then click **OK**.

Test Services is set up to run under Local System account and does not require interaction with the desktop.

NOTE

Agilent Test Services service depends on the Agilent OpenLab Shared Services (OLSS) service, as well as most of the services of OpenLab CDS software to perform operations. It will show registered service dependency on the OLSS service. In cases of system reconfiguration, make sure that the OLSS service and Agilent Test Services service are restarted to utilize the new configuration.

Licensing

Users do not need any license to log in to Test Services, execute the OpenLab Software Installation Verification Test (SVT), System Report, Connectivity test, view last test results, or access reports from previous test executions. However, to execute the OpenLab Security Test and/or the OpenLab Storage System Test, the QlaSecurity license must be available. To execute the Workflow test, the QlaCDSWorkflow license must be available. Each machine where a licensed test is executed requires one license.

The license is acquired at the first execution of a licensed test and remains acquired after the test execution completes. (It is NOT released to the license pool after the test completes.)

QualA licenses are installed from the Licenses option available in the OpenLab Control Panel Administration tab. You must install the license before you run a licensed test. If you find that the number of licenses you have is insufficient, contact Agilent support to obtain extra licenses.

Test Services utilizes the same subsystem for licensing as OpenLab CDS. In the license file, the feature name for the Security test and Storage System test is QlaSecurity and for the Workflow test the name is QlaCDSWorkflow.

NOTE

The OpenLab CDS 60-day startup license is not sufficient to run the OpenLab licensed tests in Test Services. Specific QualA licenses must be purchased and installed for licensed tests to work.

If you want to retire a server, uninstall Test Services for the QualA license to be returned to the pool of available licenses.

For tests that require a QualA license, the license is acquired when the user clicks the Start button for a licensed test. A "License Acquired" message appears next to the test name.

If there are not enough licenses available for the test execution, the test will fail at the second step "Checking for license <license name>."

You will still be able to view previous Test Services results.

Setting User Account to Run Test Services

Test Services uses OpenLab Shared Services (OLSS) authentication. For details on how to configure different authentication providers, see the *OpenLab CDS Client and AIC Guide* or the *OpenLab CDS Workstation Plus (With Content Management) Guide*.

NOTE

For Workstations without authentication, no user login is required, and this section does not apply.

Test Services user roles and privileges

After installation, log in as an OLSS Admin user. At first login, new Project roles Test Services Administrator and Test Services User are created. New privileges, Run Test Services and Manage Test Services, are created in OLSS. Test Services Administrator has all privileges from Test Services User role plus the Manage Test Services privilege. By default, any user with Everything role will have these privileges assigned automatically.

Users who do not have the Everything role will need to have the Test Services User role added to their profile through the OpenLab CDS Control Panel. At a minimum, the Test Services User role allows users to log in to Test Services and execute the System Report and OpenLab Software Installation Verification Test (SVT).

For information on adding users or updating user roles and privileges, see the OpenLab CDS Control Panel help.

For Test Services privileges using OpenLab ECM, see **[“Test Services user role creation in OpenLab ECM 3.x”](#)** on page 17.

Table 1 Privileges included with Test Services roles

Privilege	Test Services Administrator	Test Services User
View project or project group	x	x
Manage project or project group	x	x
Edit content of project	x	x
Run test services	x	x
Manage test services	x	

Required roles and privileges

In addition to the Test Services User Role, the following privileges are required to run the tests.

NOTE

For Workstation with *Authentication: None*, no privileges apply.

Table 2 Privileges required to perform tests (in addition to Test Services User Role)

Privilege	Software Installation Verification Test	System Report	Security Test	Storage Test	Connectivity Test	Workflow Test
Manage security			x	x*		
View Activity Log			x	x [†]		x
View instrument or location					x	x
Manage instrument or location					x	x
Create and modify sequence						x
Run instrument						x

Table 2 Privileges required to perform tests (in addition to Test Services User Role)

Privilege	Software Installation Verification Test	System Report	Security Test	Storage Test	Connectivity Test	Workflow Test
Delete Content of Project						x**
Edit users own running sequences Or Edit any users running sequences						x

* For Workstation with Internal/Domain authentication, the Manage security privilege is not required for the Storage test.

** For Workstation with Internal/Domain authentication, the Delete Content of Project privilege is not required for the Workflow test.

‡ For Workstation with Internal/Domain authentication, the View Activity Log privilege is not required for the Storage test.

Test Services user role creation in OpenLab ECM 3.x

At first log-in, the Test Services User role is created in OpenLab ECM 3.x, containing the following privileges:

- Content: File (View, Add)
- Content: Folder (View, Add)
- Content: File Filtering (Edit)

To use Test Services, users must be present (created/imported) in the OpenLab ECM 3.x and OLSS with their privileges updated to include the Test Services User role.

With these privileges, the user will be able to log in to Test Services, view history, download reports, execute SVT and System Report.

For information on adding users or updating user roles and privileges in OpenLab ECM, see the OpenLab ECM online help.

Client/server configuration with OpenLab ECM 3.x

To run the OpenLab Security Test, the user needs to also have the following privileges:

- System: Service Administration (Run)
- System: Audit Trail (View)
- System: Roles Configuration (View, Edit)
- System: Users/Groups Configuration (View, Edit)
- Manage Security (OLSS privilege)
- View Activity Log (OLSS privilege)

To run the OpenLab Storage System Test, the user needs to also have the following privileges:

- Content: File (Delete)
- Content: File Association (View, Add)
- Content: File Revisions (View)
- Content: Folder (Edit, Delete)
- Content: Folder Access Properties (Edit)
- Content: Move File (Run)
- Content: Move Folder (Run)
- Content: Rekey File (Run)
- System: Audit Trail (View)
- System: Filtering Configuration (Edit)
- System: Roles Configuration (View, Edit)
- System: Users/Group Configuration (View, Edit)
- System: Quick Search (Run)

To run the Workflow Test, the user needs to also have the following privileges:

- Content: File (Delete)

To run the Connectivity Test, the user needs to also have the following privileges:

- View instrument or location (OLSS privilege)
- Manage instrument or location (OLSS privilege)

Logs

Viewing the Test Services logs

The Test Services log files are at `C:\ProgramData\Agilent\LogFiles`. Note that ProgramData is a hidden folder.

The TestServices.log file contains information regarding operations performed by Test Services and its plug-ins. For easier troubleshooting, each plug-in logs the events with tags indicating the test that generated the events.

Another log file can be found on the machine where Test Services Central Management is installed, at: `C:\ProgramData\Agilent\LogFiles\TestServices_cm{timestamp}.log`. This log can be used to troubleshoot the Test Services central management services.

NOTE

When upgrading from a version of QualA prior to v3.4, path names are changed from QualA to Test Services. When upgrading using a custom path, a Test Services folder is created in the custom path.

Configuring Test Services

This section describes configuration actions an administrator can perform.
After every change, restart the Agilent Test Services service to apply the changes.

Modifying the port number

A configuration tool is available that lets you change the port for Test Services.

- 1 Go to `C:\Program Files (x86)\Agilent Technologies\Test Services`.
- 2 Run the `Agilent.TestServices.ConfigTool.exe` as an administrator.
The currently configured port is displayed along with the status of the Test Services service.
- 3 In **Test Services TCP Port**, type the new port number.

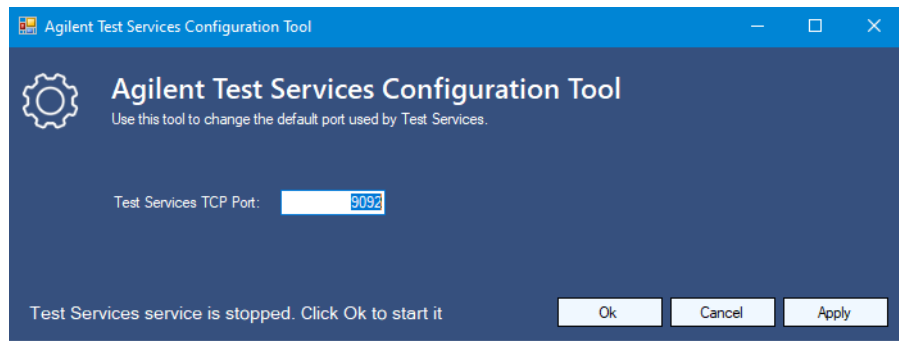


Figure 5. Test Services Configuration Tool

- 4 Click **Apply**.
The Test Services service is stopped and then started again.
- 5 Click **Cancel** to close the configuration tool.

- 6 For Client: Log in to Test Services using the new port.
(https://localhost:<new_port>). For other machines, log in to Test Services as usual (https://localhost/testservices.)
- 7 When completed, the change is logged into the following log file: C:\ProgramData\Agilent\InstallLogs\TestServicesConfigToolLog.log.
- 8 If the port is changed on a system where the Reverse Proxy is installed and running (that is, everywhere except on the Client) the reverse proxy configuration must be modified if the port has changed. The procedure is as follows.

- a As a computer administrator, open "%ProgramFiles(x86)%\Agilent Technologies\OpenLab Reverse Proxy\Apache24\conf\httpd.conf" in a text editor (for example, notepad) and modify the following section:

```
#<testserviceswebui>

ProxyPass /testservices https://localhost:9092/testservices
retry=3 acquire=3000 keepalive=On

ProxyPassReverse /testservices
https://localhost:9092/testservices

#</testserviceswebui>

#<testservicesrestapi>

ProxyPass /openlab/ca https://localhost:9092/openlab/ca
retry=3 acquire=3000 keepalive=On

ProxyPassReverse /openlab/ca
https://localhost:9092/openlab/ca

ProxyPass /openlab/testservices
https://localhost:9092/openlab/testservices retry=3
acquire=3000 keepalive=On

ProxyPassReverse /openlab/testservices
https://localhost:9092/openlab/testservices

#</testservicesrestapi>
```

Change the default port number 9092 in all 6 places in this configuration file (to the port number configured in previous steps with the tool).

- b Save the file.
- c Restart the Apache2.4 service.

NOTE

If reconfiguring a client system to port "XXX", use https://localhost:XXX/ to access test services locally. To access test services remotely, use https://<client-fqdn>:XXX/. If there is a firewall on the client system, make sure that port XXX is open.

If OpenLab CDS configuration changes after installation of Test Services

If certain changes are made to OpenLab CDS configuration after Test Services is installed, it is necessary to run a tool that creates the Test Services User role and Run Test Services privilege in the system:

```
C:\Program Files (x86)\Agilent Technologies\TestServices\
Agilent.TestServices.PrivilegeCreatorTool.exe <username>
<password> <domain>
```

The following configuration changes require running the tool:

- After installing Test Services on a CDS AIC/Client, the CDS AIC/Client is reconfigured to be connected to another OLSS server.
- After installing Test Services on a Workstation with Authentication set to **None**, the Authentication provider is changed to **Internal** or **Domain**.
- After installing Test Services on a client/server system using OpenLab ECM 3.x with OLSS Server configured to use one ECM account, the OLSS Server is reconfigured to use another account in ECM.

If authentication changes after installation of Test Services

After a change is made to authentication, you will need to log into Test Services as an admin user. This automatically updates current schedules, notifications, and parameters.

Location of Test Services Test Results

OpenLab Server and ECM XT

For OpenLab Server and ECM XT, the results of Test Services tests are saved in Content Management, in the project folder created by Test Services: **Content > Test Services Project > Test Services Results**. Each folder within the Test Services Results folder contains the files for a given test execution.

NOTE

If you upgraded from a previous version of QualA, your previous QualA result folders will remain. Tests run after the upgrade will be located in **Content > QualA Project > QualA Results**.

ECM 3.5 and 3.6

For ECM 3.5 and 3.6, the results of Test Services tests are saved in ECM, in the project folder created by Test Services: **Content > TS > TS > Test Services Project > Test Services Results**. Each folder within the Test Services Results folder contains the files for a given test execution.

Security and Certificates

By default, Test Services uses secure communications (HTTPS) on any machine where Test Services is installed.

There are some situations where installation of a third party certificate is necessary for access to Test Services over a secure connection:

- On devices where Test Services is not installed but secure access to the Test Services URL on another machine is required.
- Situations where use of custom or third party certificate is required.

If you access the Test Services URL from a device that does not have a valid certificate, you will see a warning that the connection is not private. You can still access Test Services without the certificate installed, however you will get another warning once you open Test Services.

Use the following procedure to install a third party certificate and then configure Test Services to use it.

Configure Test Services to use a third party certificate

If you want to use an installed third party certificate, you must configure Test Services to use it.

NOTE

Before using this procedure, obtain and install your third party certificate.

- 1 Locate and copy the serial number of the certificate you want to use.
 - a From the Windows Control Panel, in Administrative Tools, go to **Manage Computer Certificates**.
 - b In the **Personal** folder, locate the certificate you want to use.
 - c Double-click the certificate and select the **Details** tab.
 - d Select **Serial number** and copy it.

Configure Test Services to use a third party certificate

- 2 As an administrator, in the Command Prompt, run the Test Services configuration tool with the -cert option, and the serial number you copied. The following is an example, where the serial number is 00ECF5187D8D286484.

```
C:\Program Files (x86)\Agilent Technologies\Test  
Services>Agilent.TestServices.TestServicesConfigTool.exe  
-cert 00ECF5187D8D286484
```

- 3 To test the secure connection, log in to Test Services:
<https://localhost/testservices> (or <https://localhost:9092> on Client machines).
If the connection is secure, the security icon to the left of the Test Services URL in your browser will appear "locked."  <https://>



3 Troubleshooting

Troubleshooting 27

Error Messages 31

This chapter contains information on how to troubleshoot issues and resolve error messages that occur while using Test Services.

Troubleshooting

Results reported as not passed

The main task for Test Services is to ensure that the OpenLab software operates as expected. It is therefore expected that the resulting report always produces the predefined values with a result of "Passed." However, in rare occasions when the software is not performing as expected, the report will show status of "Failed" or "Not Available," in which case an administrator can perform corrective actions.

Failed report status

Test Services can show a "Failed" report status for various reasons, such as (but not limited to):

- Corrupted OpenLab CDS files
- Windows Update was installed that is not compatible with OpenLab CDS

To mitigate these issues, perform the following actions:

- To verify the software (both OpenLab CDS and Test Services) was not modified incorrectly, run the software verification tool that is preinstalled with OpenLab CDS.
- Run OpenLab CDS on the machine. Launch the OpenLab Control Panel, launch an instrument, and try to perform a run. This should highlight whether there are any issues with CDS.
- On rare occasions, a Windows Update may introduce non-backward compatibility of the software that is based on .NET Framework. If you notice that this is the case, try to identify the problematic Windows Update and rollback its changes.

CAUTION

Uninstalling a Windows update may have security or other implications. Before you uninstall the Windows update, make sure that you understand the implications and the effect that uninstalling might have on the system. Perform a system restore checkpoint and verify that you have options to restore the system to the state prior to uninstalling an update.

Connectivity Test failed

"Failed" status of the Connectivity test appears when at least one component is not connected.

- If executed against an AIC, "Failed" status of the Connectivity test appears when at least one instrument is not connected.
- If executed against a Client, "Failed" status of the Connectivity test appears when at least one AIC in the system is not pingable from the Client.

Report status Not Available

This status is more common than the "Failed" result. It does not necessarily mean that something is wrong with OpenLab CDS, but if it occurs regularly it should be verified that the operations of OpenLab CDS are not impaired.

"Not Available" result is shown in cases where:

- The test was not run to its completion (due to service restarts or similar.)

Sequence hangs during workflow test

When running the Workflow test where the machine is in a domain, the Workflow test hangs at step "Starting sequence <sequence>."

To execute the Workflow test successfully, the Test Services service must run under the same user account as Instrument service.

NOTE

The user may or may not be a domain account. In case of a workstation that is not in a domain, this user may just be a local windows user account.

Perform the following steps to update the Test Services service to use the same user that is used for the Instrument service:

- 1 Verify the user account under which the Instrument service is running.
- 2 Log in as Windows domain user who has local administrative privileges.
- 3 Go to **Control Panel > All Control Panel Items > Administrative Tools** and double-click **Services**.
- 4 Right-click **Agilent Test Services Service** and select **Properties**.
- 5 Click the **Log On** tab, select **This account** and enter the login credentials of the user that is used for the Instrument Service. Click **OK**.

- 6 The Services window confirms that the account has been granted the **Log On As A Service** right. Click **OK**.
- 7 Click **OK** in the Services window to acknowledge that **The new logon name will not take effect until you stop and restart the service.**
- 8 Restart the Agilent Test Services Service.
- 9 Log in using the same user account specified in the **Log On** tab of Agilent Test Services Service Properties window.

Test Services projects and instruments

The Test Services Project is created during the first execution of any test. During the Workflow test, the Test Services project is configured with additional parameters, and the required Test Services instrument is created. If for any reason these are deleted or modified, Test Services will recreate them on the next run.

Email notification not working

If email notifications are not working, use the following procedures to troubleshoot.

- Check to make sure that email is set up correctly in OpenLab CDS Control Panel. If it is not set up, the email notifications will fail. For information on how to set up email, see the OpenLab CDS Control Panel online help.
- Send a test email from OpenLab CDS Control Panel to make sure that your email settings are working.
- Test the email from Test Services.

a Set up email notifications in Test Services.

b In the browser, type the following URL. For client/server systems, the server is the address of OpenLab Server, or an OLSS server with ECM 3.x backend. For Workstation or Workstation Plus, this is the machine itself:

```
https://<server:port>/openlab/testservicesserver/v1/health/email
```

For example,

```
https://localhost:52088/openlab/testservicesserver/v1/  
health/email
```

This sends an email using default values from the Notification page in Test Services. It uses the SMTP settings from OpenLab Control Panel.

A test email is sent with the following subject: Test Services Test email.

NOTE

Email can only be sent from server, Workstations or Workstation Plus machines (no clients or AICs.)

Error Messages

The following lists all error messages that Test Services can generate, with actions that can be taken to mitigate the issues.

Test Services error messages

Account is expired

- Verify that the user account is valid in the OpenLab CDS Control Panel.

Test Services could not acquire a license. Please contact your system administrator to resolve the issue.

- Verify that you have obtained a valid license for QualA in OpenLab CDS Control Panel, and that the license hasn't expired.
- Verify that you have enough licenses available. If you find that number of licenses is insufficient, contact Agilent support for additional volumes.
- Verify that the Flexera software (third party system for managing licenses) is operational by opening its dashboard. (By default, it can be found on <http://localhost:8090> for Server or <http://localhost:8095> for Workstation and WorkstationPlus, where localhost denotes a machine with the licensing information.) You should see a feature named QlaSecurity, with quantity matching the number of licenses obtained.
- Restart the following Agilent OpenLab Windows services: License Server, Licensing Support, Shared Services, Test Services. (To open the services page, press Windows button + R, type `services.msc` and then press **Enter**.)

Test Services Service is down or unavailable - please contact system administrator for assistance

- This error occurs when the Test Services windows service stops being reachable – either because of network errors, or the service itself stopped. Make sure that the port is still open, and that the service is up and running. If necessary, restart the Windows Service (Agilent Test Services, in the `services.msc`).

Bad username or password

- This is an exception thrown by the underlying authentication mechanism of the Shared Services. Make sure that the user has correct user name and password in the OpenLab CDS Control Panel administration page.

Logon credentials are either invalid or expired

- Troubleshooting steps are the same as for [Bad username or password](#).

Authentication provider is not available

- This is an exception thrown by the underlying authentication mechanism of the Shared Services. Make sure that the OpenLab CDS Control Panel does not report any errors and resolve the issues there.

Connection to server has failed

- Troubleshooting steps are the same as for [Authentication provider is not available](#).

Failed to validate credentials

- Troubleshooting steps are the same as for [Authentication provider is not available](#).

OLSS service is not available

Troubleshooting steps are the same as for [Authentication provider is not available](#).

Test Services cannot continue verification - please contact system administrator for assistance

- Troubleshooting steps are the same as for [Authentication provider is not available](#).

User has empty password

- Troubleshooting steps are the same as for [Authentication provider is not available](#).

User is disabled

- Troubleshooting steps are the same as for **Authentication provider is not available**.

User is temporarily blocked

Troubleshooting steps are the same as for **Authentication provider is not available**.

User must change password

Troubleshooting steps are the same as for **Authentication provider is not available**.

User is not mapped but is authenticated

- Troubleshooting steps are the same as for **Authentication provider is not available**.

Insufficient privileges to run Test Services

- User that logged in does not have a privilege to access the Test Services. Make sure that the user has the role Test Services User. (See **“Setting User Account to Run Test Services”** on page 15.)

Test Services cannot create Test Services Project - please contact system administrator for assistance

- This error can occur if the content management is unavailable, or it is available but there are some other issues in the content management (for example, disk space, licensing issues, or similar). Make sure that there are no issues with the content management. Also, you can try to remove the Test Services project from the Control Panel and it will be created automatically on the next run of Test Services. All previous results will remain in the Content management untouched.

Report is not available in the storage

- The error shown means that the PDF report was not found in the content management designated folder. Make sure that the current user has permissions to the folder specified in the configuration of Test Services. This message can also occur if the file was removed or archived directly using the content management. In this case, restart the Test Services execution.

Email server settings are not configured in Control Panel for OpenLab software

- Test Services uses email settings from OpenLab Shared Services to send email notifications. If you see this error, go to the OpenLab Control Panel and set up your email preferences.

OpenLab CDS Workflow Test error messages

Error configuring Test Services instrument - please contact system administrator for assistance

- This error can occur if the instrument is currently in use. Make sure it is not currently in use. If this fails, delete the instrument, as it will be created automatically on the next run of Test Services.

Test Services cannot create instrument - please contact system administrator for assistance

- Troubleshooting steps are the same as for [Error configuring Test Services instrument - please contact system administrator for assistance](#).

Error configuring Test Services Project - please contact system administrator for assistance

- This error can occur if the Content Management is unavailable, or it is available but there are some other issues in the Content Management (for example, disk space, licensing issues, or similar). Make sure that there are no issues with the Content Management. Additionally, you can try to remove the project from the Control Panel and it will be created automatically on the next run of Test Services. All previous results will remain in the Content management untouched.

Test Services project is required for Test Services verification - please contact system administrator for assistance

- This error occurs if the instrument is found, but the project is not. Since Test Services creates both automatically, this message only occurs if somebody interfered with the process of Test Services during the execution of the steps. Make sure that Test Services is uninterrupted from the OpenLAB CDS, and then retry Test Services.

Valid local instrument controller is required for Test Services verification - please contact system administrator for assistance

- This occurs if the Instrument Controller is unavailable while configuring the Test Services instrument. Verify that the AIC is operational and retry.

Error submitting sequence - please contact system administrator for assistance

- Sequence run can fail for any number of reasons. Refer to the Test Services logs found in the OpenLAB CDS Control Panel > Administration > Diagnostics section for further information.

In This Book

Administration for Agilent Test Services

- QualA Licensing
- User account set up
- Troubleshooting

www.agilent.com

© Agilent Technologies, Inc. 2022
Document No: D0015137
April 2022

