



Test Services

User Guide

Notices

Document Identification

DocNo D0146323 Rev. A.00
04/2026

Copyright

© Agilent Technologies, Inc. 2022-2026

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This guide is valid for the 3.7 revision or higher of the Test Services program until superseded.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 document provides instructions on how to use the Test Services program.

Test Services is installed automatically when you install OpenLab CDS. It includes a suite of verification tests that are available without a license. Test Services for OpenLab CDS provides licenses that enable additional CDS verification tests.

1 Introduction

This chapter provides an overview of how Test Services works.

2 Using Test Services

This chapter provides information on how to log in, run verification tests, and view reports. It also describes how to schedule tests and set up credentials and email notifications.

Table of Contents

1	Introduction	5
	Overview	6
2	Using Test Services	8
	How to Use Test Services	9
	Licensing for Test Services tests	9
	Logging in to Test Services	9
	Test Services navigation	12
	Running Test Services Tests	13
	OpenLab Software Installation Verification Test	15
	System Report	16
	OpenLab Security Test (License Required)	16
	OpenLab Storage System Test (License Required)	18
	OpenLab CDS Workflow Test (License Required)	19
	OpenLab CDS Connectivity Test	20
	Reviewing Test Services Reports	22
	Test report contents	25
	Aborting Test Execution	28
	Scheduling Test Services Tests	29
	Create a test schedule	29
	Enable or disable a test schedule	33
	Change a test schedule	33
	Delete a test schedule	34
	Settings	35
	Setting up credentials for running tests	35
	Setting up email notifications	36



1

Introduction

Overview 6

This chapter provides an overview of how Test Services works.

Overview

Test Services provides a simple and fully automated mechanism for verification of software and system functionality. Test Services consists of the Test Services framework, with optional plug-ins that add verification tests. When run, Test Services provides a series of user-selectable tests that confirm the system is operating as intended. Test Services tests are listed in [Table 1](#). Some tests are installed and enabled by default without a license. Some tests are installed, but require a license in order to run them.

NOTE

When you run Test Services on a given system component (AIC, client, server) only tests that are available on that component are accessible. If a test is part of a plug-in that is not installed, it does not appear.

Table 1. Available tests in Test Services

Test	Tasks performed	Workstation	Workstation Plus	Server/ECM XT	Client	AIC
System Report	• Lists installed applications and patches • Indicates changes since the last system report was executed.	Available	Available	Available	Available	Available
OpenLab Security Test	• Confirms password policy enforcement • Confirms roles-based access enforcement • Verifies activity logs	Available (Internal/Domain authentication) or Disabled - no authentication) (license required)	Available (license required and included)	Available (license required)	Not recommended	Not recommended
OpenLab Storage System Test	• Verifies storage system's functionality for storing, securing, and maintaining files and folders	Available (license required)	Available (license required and included)	Available for OpenLab ECM and ECM Server/XT (license required)	Not recommended	Not recommended

Table 1. Available tests in Test Services (continued)

Test	Tasks performed	Workstation	Workstation Plus	Server/ECM XT	Client	AIC
OpenLab Software Installation Verification Test	Verifies integrity of installed Agilent Software binaries by running the Software Verification Tool (SVT)	Available	Available	Available	Available	Available
Workflow Test (Requires OpenLab CDS Plug-in)	Creates a project and virtual (Data Player) instrument, submits a sequence and verifies the software algorithms	Available (license required)	Available (license required and included)	Not available	Not available	Available (license required)
Connectivity Test (Requires OpenLab CDS Plug-in)	- Network connectivity to all system AICs - Connectivity to instruments that are configured on the AIC	Available	Available	Not available	Available	Available

It is also possible to create a test schedule for automatic execution of selected tests. If applicable, you can also select where to run the tests (for example, on a client or AIC.)

You can also set up notifications to send emails to selected users to notify them of test execution and results.

2 Using Test Services

How to Use Test Services 9

Licensing for Test Services tests 9

Logging in to Test Services 9

Test Services navigation 12

Running Test Services Tests 13

OpenLab Software Installation Verification Test 15

System Report 16

OpenLab Security Test (License Required) 16

OpenLab Storage System Test (License Required) 18

OpenLab CDS Workflow Test (License Required) 19

OpenLab CDS Connectivity Test 20

Reviewing Test Services Reports 22

Test report contents 25

Aborting Test Execution 28

Scheduling Test Services Tests 29

Create a test schedule 29

Enable or disable a test schedule 33

Change a test schedule 33

Delete a test schedule 34

Settings 35

Setting up credentials for running tests 35

Setting up email notifications 36

This chapter provides information on how to log in, run verification tests, and view reports. It also describes how to schedule tests and set up credentials and email notifications.

How to Use Test Services

Test Services is started by opening an Internet browser URL for Test Services installed on an OpenLab CDS AIC, Client, Server, Workstation, or Workstation Plus.

For server, AIC, Workstation Plus, and workstations, use

`https://hostname.domain.com/testservices`

or

`https://localhost/testservices`

On client systems, use

`https://localhost:52088/testservices` to access test services locally.

To access test services remotely, use

`https://<<client-fqdn>>:52088/testservices`

If there is a firewall on the client system, make sure that port 52088 is open.

Licensing for Test Services tests

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. For more information, see the *Test Services Administration Guide*.

Logging in to Test Services

After you open the Test Services URL in your browser, the Test Services Log in page appears.

NOTE

If you are using a file-based Workstation with authorization set to “None,” no login is required.

Type a valid user name and password that you use to log in to OpenLab CDS (and if applicable, a Domain), and then click **Sign in**.

NOTE

For the first login after installation, the user should have OpenLab Shared Services administrator role, to allow the framework to finish registering itself with the OpenLab Server.

After the first login, the Test Services User role is sufficient to log in, execute the Software Installation Verification Test and System Report, and view previous results in history.

To execute the OpenLab Security Test and OpenLab Storage System Test, the System Administrator role with associated Manage Security privilege is required.

For information on editing a user account to add privileges, see the OpenLab CDS online help and the *Test Services Administration Guide*.

After you log in, the Test Services home page appears. Tests that do not require additional license are displayed at the top. Tests that require a license are displayed under **Licensed Tests**.

By default, only recommended tests are displayed. To display all available tests, whether or not they are recommended on the machine from which you run Test Services, click **Show all tests**. See “Available tests in Test Services” on page 6.

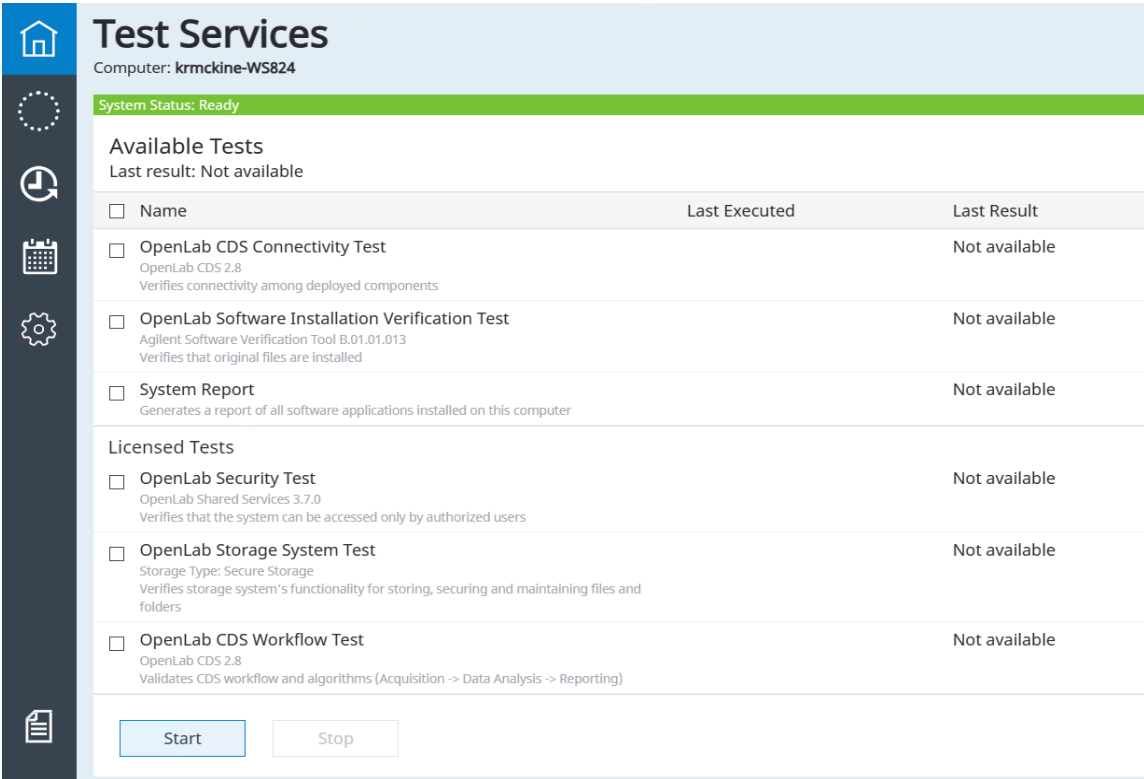


Figure 1. Test Services home page example

NOTE

Where a license is required, it must be installed before running the test. The license is acquired at the first execution of the test, and the display shows License Acquired.

The license remains acquired after the Security or OpenLab Storage System Test execution completes. (It is NOT released to the license pool after the Security or OpenLab Storage System Test completes.)

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

Test Services navigation

On the left side of the home page is a navigation pane where you can select the following task areas:



Home:

Displays list of user-selectable tests. After selecting desired tests, you start the tests from this window. You can also view last execution results and access reports of the last execution.



Execution Details

Displays the current test execution in progress with detailed progress steps and their status.



Execution History

Displays list of tests that were performed. You can select a date range or create a filter and open reports.



Test Schedules

Displays a list of current test schedules. You can create, edit, enable, disable, or delete test schedules.



Settings

Displays configurable settings for Test Services

Credential Settings - you can configure credentials for running ad-hoc and scheduled tests.

Notification Settings - you can enable, disable, and configure notifications for test executions. By default, notifications are disabled.



Help and documentation

Links to Test Services User Guide and Administration Guide.



User log out. User is automatically logged out due to inactivity after the timeout defined in the OpenLab Shared Services Security Policy. Note: For file-based workstations with authentication set to "None," this button is not displayed.

Running Test Services Tests

NOTE

Test Services only enables the tests when the minimum prerequisites for the test are met. For example, the Software Installation Verification Test is disabled if the Software Verification Tool (installed with OpenLab CDS) is not present on the system.

If plug-ins are installed, those tests appear in **Available Tests**.

To start verification tests,

- 1 In the home page, select the boxes for the tests you want to perform (**Figure 2**). To select all tests, click the box next to **Name**.

NOTE

You can select tests and start their execution when the System Status is *Ready*. If *Running* or *Aborting*, you must wait until the execution finishes.

- 2 Click **Start**.

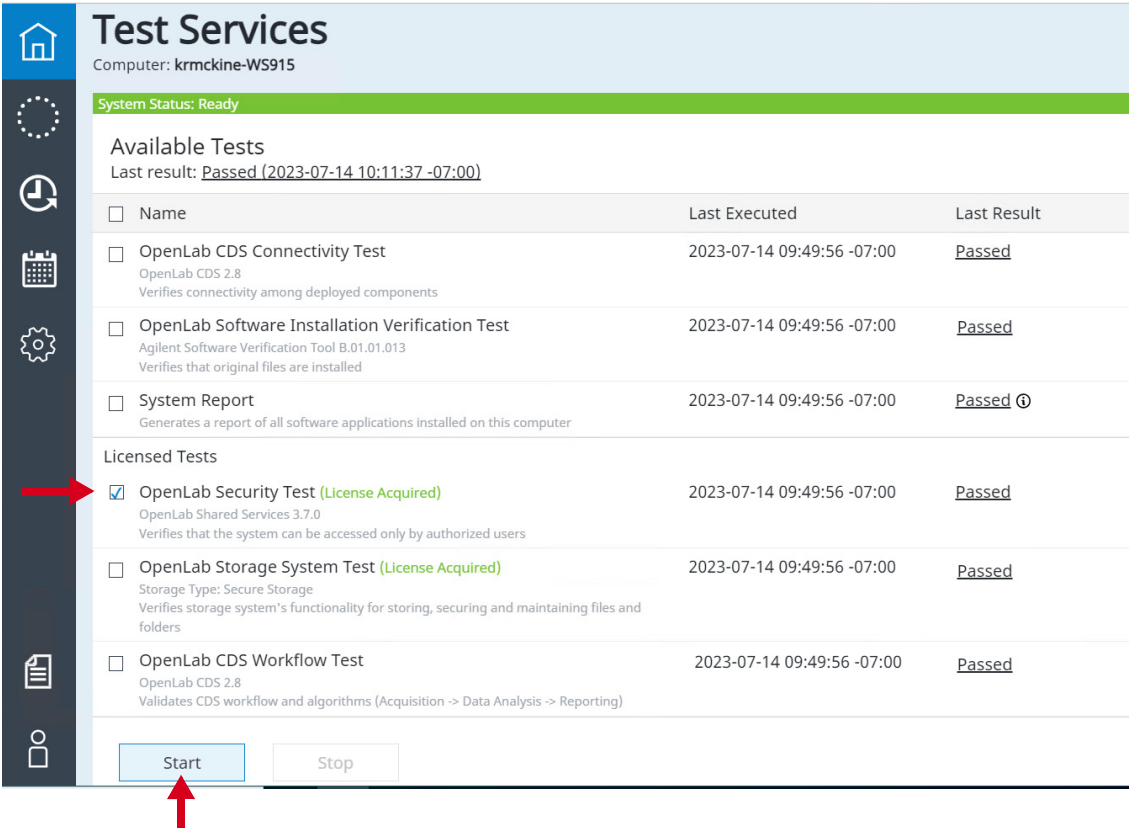


Figure 2. Start tests from the home page

When a test starts, the system status at the top of the home page changes to *Running*, along with additional information such as the number of tests queued for execution.

NOTE

A "Passed with additional information" icon [Passed](#) ⓘ appears in cases where a test passed, but there is additional information to review in the report.

3 To view the status of the selected tests, select the **Execution Details** page.

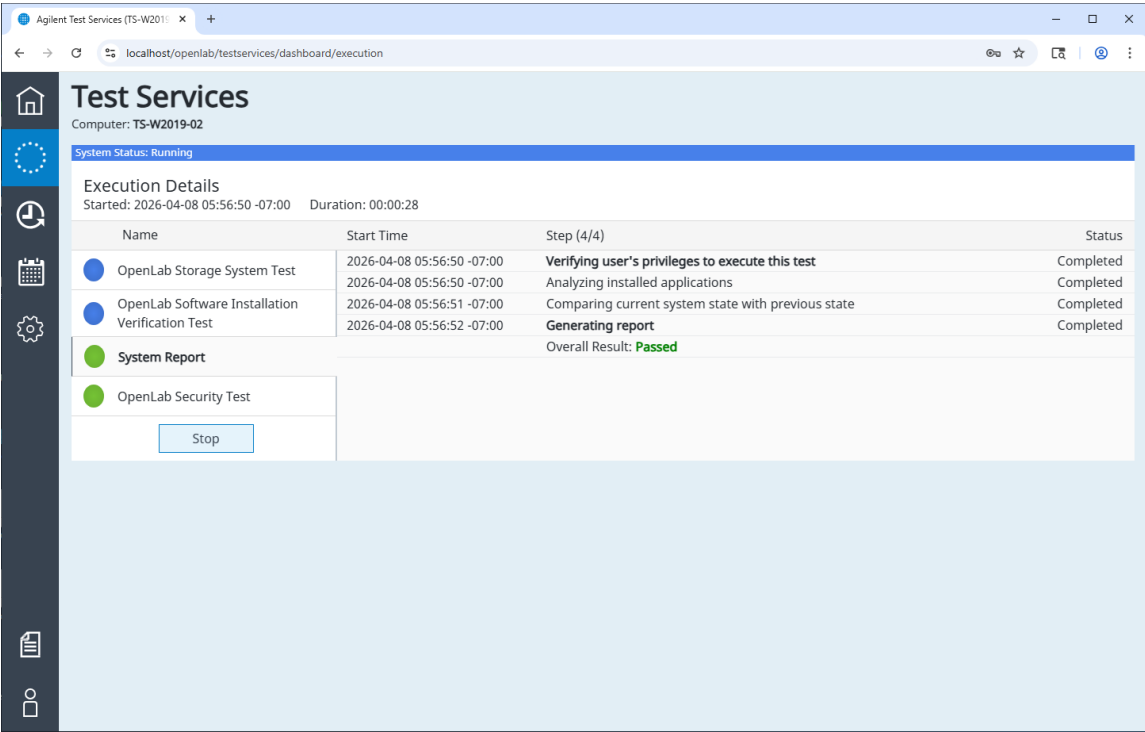


Figure 3. Execution Details shows progress of tests

- 4 To see the progress/status of a test (Pending/In Progress/Completed/Aborting) click the test.
- 5 To abort the tests, click **Stop**.

A description for each test is given in the following section.

OpenLab Software Installation Verification Test

Verifies the integrity of installed Agilent software binaries by running the Software Verification Tool (SVT). This test is available for all systems.

System Report

Generates a System Report including all software applications installed on this computer.

- System Details - hardware and operating system configuration, including antivirus installed, time synchronization information, and information on users who are allowed to change date/time and time zones
- Software Applications - installed software applications including version, publisher, and date installed
- Updates - installed updates including type and date installed
- Highlights changes from the previous time the System Report was generated

If a change was detected since the last system report, a “Passed with additional information” icon appears. **Passed** ⓘ

Warning messages in the report are generated for critical or unexpected changes (for example, antivirus is not up to date or installed, if Windows “Users” group can change date/time or time zone, or more than 24 hours elapsed since time was synchronized.)

This test is available for all systems.

NOTE

Antivirus information is not available on Windows Server systems.

OpenLab Security Test (License Required)

The OpenLab Security Test verifies the system can only be accessed by authorized users.

NOTE

If you are using Domain authentication provider, you must configure the Security Test before it will be available. For more information, see **“Configuring the OpenLab Security Test”** on page 17.

For this test, two users are created and imported to the system: admin and non-admin.

Then, using the created/imported users, the security tests are executed in OpenLab Shared Services (OLSS).

Configuring the OpenLab Security Test

If you are using Domain authentication provider, the OpenLab Security Test needs to be configured before it becomes available for execution. In this case, you will see a **Configure** button below the OpenLab Security Test in the home page. Click the button to open a Test Configuration form. In the form, enter user credentials for test users (admin and non-admin user) that the test will use for the test execution.

You can also select to remove the test users from the system at the end of the test. The entered data can also be saved for later use.

CAUTION

If "Remove users at the end of the test" is selected, and the Security and Storage tests are run simultaneously, the Storage test might fail with the message "Account is not authorized in this system. Contact system administrator." This is caused when the users from the Security test are deleted before the Storage test is completed. This can also happen if the test is run from two different machines at the same time. To avoid this, set different users for the Security and Storage tests.

Configuration is performed once, and applied globally to all machines in a client/server environment.

NOTE

If a currently logged-in user is set as a test user, and you select to remove the test users at the end of the test, the test will fail. The system does not allow removal of test users currently logged into the system. (For example, removing yourself from the system creates an error.)

NOTE

In client/server environments, when the Test Services administrator sets this information on one machine, it will appear on all machines within 3 minutes.

OpenLab Storage System Test (License Required)

This test verifies functionality for storing, securing, and maintaining files and folders. The test performs slightly different verifications depending on what type of storage system you are using.

NOTE

If you are using Domain authentication provider, you must configure the OpenLab Storage System test before it will be available. For more information, see [“Configuring the OpenLab Storage System Test”](#) on page 18.

Configuring the OpenLab Storage System Test

For Secure Storage, if you are using Domain authentication provider, the OpenLab Storage System Test needs to be configured before it becomes available for execution. In this case, you will see a Configure button below the OpenLab Storage System Test in the home page.

Configuration is performed once, and applied globally to all machines in a client/server environment.

For Secure Storage with Domain authentication

- 1 Click **Configure** to open the Test Configuration form.
- 2 Enter user credentials for test users (Test Services admin and non-admin user) that the test will use for test execution.
- 3 Select to remove the test users from the system at the end of the test. (Optional)
- 4 Click **OK**.

NOTE

In client/server environments, when the Test Services administrator sets this information on one machine, it will appear on all machines within 3 minutes.

For OpenLab ECM as a storage provider

- 1 Click **Configure** to open the Test Configuration form.
- 2 By default, **Verify quick search by uploading a file and searching for it** is selected. The test must wait for the file to be processed and indexed before the test can search for it. You can change the time to wait for the processing and indexing of the file before timing out. (Default is 300 seconds, maximum value is 9999 seconds)
- 3 Click **OK**.

OpenLab CDS Workflow Test (License Required)

During the OpenLab CDS Workflow Test, Test Services creates a project and virtual (data player) instrument and submits a sequence to verify the software algorithms.

Test Services uses sets of reference files, which are valid for the current version of OpenLab CDS. All necessary files, including report templates, sequences, acquisition and processing methods, are deployed together when Test Services is installed.

At the end of the workflow tests, a Test Services Algorithm Evaluation and OpenLab CDS Workflow Test Report are created. A summary of the Test Services Algorithm Evaluation report is also included in the Summary Report.

License required for OpenLab CDS Workflow Test

To execute the Workflow test, the QlaCDSWorkflow license must be available. Each machine where the Workflow test is executed requires one QlaCDSWorkflow license. The license is acquired at the first execution of the Workflow Test and remains acquired after the Workflow Test execution completes. (It is NOT released to the license pool after the Workflow Test completes.)

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

Using Test Services

OpenLab CDS Connectivity Test

Test Services licenses are installed from the Licenses option available in the OpenLab Control Panel Administration tab. If you find that the number of licenses you have is insufficient, contact Agilent support to obtain additional licenses. For details on how to install OpenLab licenses, see the OpenLab Control Panel online help.

Required roles and privileges

For information on the roles and privileges required to run this test, see the Required roles and privileges topic in the *Test Services Administration Guide*.

OpenLab CDS Connectivity Test

On a client, this test checks connectivity to the AICs in the system. On an AIC, Workstation, and Workstation Plus, this test checks for connectivity to instruments that are configured on the AIC.

Required roles and privileges

For information on the roles and privileges required to run this test, see the Required roles and privileges topic in the *Test Services Administration Guide*.

Connectivity tasks performed

The OpenLab CDS Connectivity Test performs the following:

- Confirms network connectivity between components: Client > AIC.
- On a client, this test checks connectivity to the AICs in the system. On an AIC, Workstation, or Workstation Plus, this test checks for connectivity to instruments that are configured on the machine.
- Not applicable to server.

NOTE

In case some instruments are not connected, the following message is displayed “One or more instruments are not connected to this AIC,” including the list of instruments not connected to the AIC.

If this message appears, connect any disconnected instruments prior to the test execution. If at least one instrument is not connected, the test will fail.

Reviewing Test Services Reports

When a test execution is completed, the test results are available in the home page or from the Execution History page, along with previous test executions.

Access reports from the home page

To view the reports from the last set of tests, click the link next to Last result: displayed under Available Tests. For example,

Last result: Passed (2018-10-05 11:39:58 +02:00)

The link opens a list of all reports generated in the last test execution, including the Summary report.

For example, if you executed in one run the Security Report and System Report, this link should show the Security Test Report, System Report, and Summary Report.

To view the reports from execution of a particular test, click the Last Result link for that test.

The link opens a list of all reports generated by this particular test (From the preceding example, clicking the **Security Test** link shows only Security Test report, no Summary Report).

Access reports from the Execution History page

The Execution History page provides a table containing test results. Included are the date and time the test was executed, the user who executed the test, the result of the test (Passed, Failed, Aborted), and the location in Secure Storage or OpenLab ECM where the reports are stored.

NOTE

For file-based workstation, the user who executed the test is not displayed.

Test Services
Computer: KM4

Execution History

Date Range
From 2019-12-31 To 2020-01-31 User Result Filter Clear

Execution Time	User	Result	Reports
2020-01-31 08:10:20 -08:00	admin	Passed	KM4-2020-07-29 01-31 08-10-20--08-00-A
2020-01-31 08:09:06 -08:00	admin	Passed	KM4-2020-07-29 01-31 08-09-06--08-00-A
2020-01-31 07:00:45 -08:00	admin	Passed	KM4-2020-07-29 01-31 07-00-45--08-00-P

Figure 4. Execution History

To filter the results

Using the displayed fields, you can filter the displayed tests using one or more of the displayed fields.

- 1 Under Date Range, select **From** and **To** dates to display only tests that were executed during that date range. By default, all statuses from all users for the last month are displayed.
- 2 Under User, type the User name who executed the test. (Not available for file-based workstations.)
- 3 Under Result, click the down arrow and select the result to apply to the filter. Choices include Passed, Failed, All, and Aborted.
- 4 To apply your selected filter, click **Filter**.
- 5 To clear the current filter criteria and set to default, click **Clear**.

To view available reports for a test

- 1 In the list of executed tests, click a test to display the available reports (Figure 5).

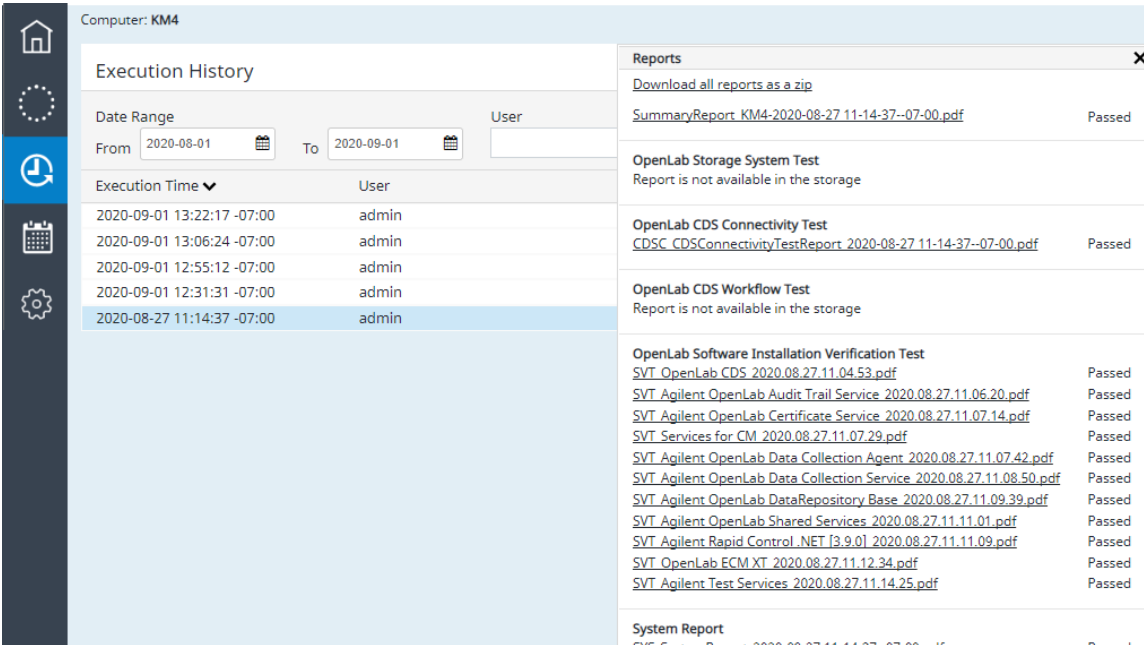


Figure 5. Available reports for selected test

The available reports are listed, along with the test result (Passed, Failed, Aborted).

- a The link at the top, under “Download all reports as a zip,” is a zip file containing all reports generated by tests in the selected test execution, including the Summary report. Click the link to download the zip file to your local machine, where you can unzip the reports and view them.
- b To view a report, click the report name. The report is downloaded to your local machine where you can open and view it.

Test report contents

The following test reports are available from the Test Services framework.

Common to all reports

- Host name, User, Date, Submitted by, Executed by (only in case of Credential users), and Overall test result (passed/failed/unavailable)

System report

This report contains version information for everything currently installed on the host machine. Any changes from the last time this report was executed are highlighted. The report contains sections for

- System Details
- Software applications
- Updates
- Summary of changes detected

OpenLab Security Test report

This report contains results of the OpenLab Security Test:

- Product name and version (if available)
- Authentication configuration details - Internal, Domain, or ECM
- Test users and roles
- Results of tests

Software Installation Verification report

For each OpenLab software component, a separate SVT report is generated that contains:

- Software application details
- Summary of results for

- File Report Summary
- GAC File Report Summary
- Files Registration Report Summary
- Registry Report Summary
- All installed files
 - Full installation path
 - Installed version

OpenLab Storage System Test report

The OpenLab Storage System Test report contains the following information:

- Time of test execution
- A user who executed the test
- Hostname
- Overall test result
- Test product and version (Local File System, OpenLab ECM, or Secure Storage)
- Product configuration (storage type, authentication provider, activity logs ON/OFF etc.)
- Test users and roles assigned
- Test configuration details
- Test procedure and steps with statuses (Passed/Failed)
- Activity logs checks

Summary report

The summary report contains a summary of the results of all the tests executed at a given date and time.

OpenLab CDS Connectivity Test report

This report is only generated when you have the OpenLab CDS Plug-in installed. It is generated when the OpenLab CDS Connectivity Test is performed.

The CDS Connectivity test report contains results of the verification of the communication between different OpenLab components and instruments.

Algorithm Evaluation report (OpenLab CDS Workflow Test)

This report is only generated when you perform the OpenLab CDS Workflow Test. It contains evaluation of all algorithms verified by the test, including expected result, observed result, and status (pass or fail for example.)

Workflow Test Report

This report is only generated when you perform the OpenLab CDS Workflow Test. It contains verifications of CDS workflow and algorithms (acquisition > data analysis > reporting.)

Aborting Test Execution

You can abort the test execution process from either the home page or the Execution Details page.

To abort the Test Services test process from the home page

- Click the **Stop** button.

To abort a test from the Execution Details page

- Click **Stop**.

The abort may not happen immediately. When it does, the current step in the table will show that the test is Aborting.

NOTE

The Stop button aborts the whole test execution—not just one test.

To stop a test, only the Run Test Services permission is required. (To start a test you need the Run Test Services permission in addition to any other permissions required by the test.)

What happens when you abort the test execution

- The System Status changes to “Aborting” during the abort process.
- The test status of Pending and Running tests change to “Aborting.” (Completed tests statuses remain “Completed.”)
- The Summary report is created with status of “Aborted.”
- The result of the tests that were not finished due to the aborted test execution shows “Not Available.”
- The result of the tests that were finished display “Passed” or “Failed” and the reports generated by these tests can be downloaded and viewed.

Scheduling Test Services Tests

You can schedule tests to run automatically on machines where Test Services is installed.

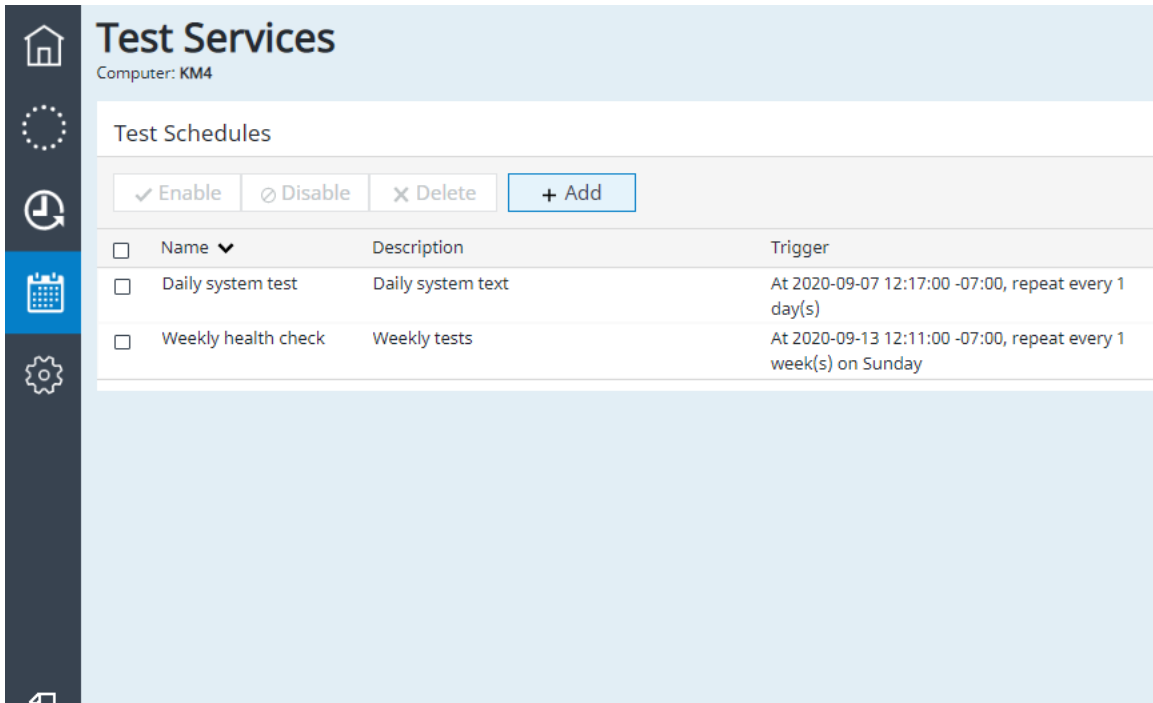
- A schedule can include one or more tests. You can include all recommended tests or a set of user-selected tests. You can also specify required parameters for the scheduled tests.
- For client/server environments, a schedule can be applied to one or more target machines, including all machines, instrument controllers, clients, servers, or specific machines.
- Tests are run using the credentials of the user who created or modified the test schedule, unless another user is specified. See “Setting up credentials for running tests” on page 35.
- Creating, editing, and deleting test schedules requires the user have the Test Services Admin role.
- Creation, modification, and deletion of test schedules are logged in the OpenLab Shared Services system log. Once a schedule is set up, the scheduled tests run whether or not you are logged into Test Services.

Create a test schedule

Test schedules are created or changed from the **Set up Test Schedule** page. This page appears when you click the calendar icon in the Test Services navigator.

If no test schedule exists, a blank test schedule appears where you can create a test schedule, using the following procedure.

If there are already one or more test schedules, a list of the test schedules appears.



The screenshot shows the 'Test Services' interface for a computer named 'KM4'. On the left is a dark sidebar with icons for Home, Status, History, Schedules (highlighted), and Settings. The main area is titled 'Test Services' and 'Test Schedules'. Below the title are buttons for 'Enable', 'Disable', 'Delete', and '+ Add'. A table lists the current test schedules:

☐	Name ▼	Description	Trigger
☐	Daily system test	Daily system text	At 2020-09-07 12:17:00 -07:00, repeat every 1 day(s)
☐	Weekly health check	Weekly tests	At 2020-09-13 12:11:00 -07:00, repeat every 1 week(s) on Sunday

Figure 6. List of test schedules

To create a test schedule, click **+ Add**, and set up the new test schedule using the following procedure.

Test Services
Computer: TS-W10-04

Set up Test Schedule

Configure one or more tests to be automatically executed.

Schedule Name:
Provide a short name for the scheduled task (<125 characters). This name will be displayed in Schedule Task tab.

Schedule Description:
Optionally, provide a detailed description for the scheduled task. This description will be displayed in Schedule Task tab.

Test Selection: ▼
Select the test(s) that need to be executed.

Machine Selection: ▼
Select the machine(s) that need to be executed.

Schedule Specify when and how often scheduled task will be executed.

[Save](#) [Cancel](#)

Figure 7. Set up Test Schedule

To set up a test schedule,

- 1 Type a schedule name.
- 2 Type a description for the schedule (optional.)
- 3 From the **Test Selection** drop-down, select what tests you want to run:
 - Run all tests** executes all available tests.
 - Run all recommended tests** executes all tests recommended for the machine where the tests are run.
 - Run specific tests** opens a menu where you can select from all tests.
- 4 From the **Machine Selection** drop-down, select from the following:
 - All machines**

Using Test Services

Create a test schedule

All Instrument Controllers

All Client Machines

Server Machine

Specific Machine(s) lets you select the machines where you want to execute the schedule.

- 5 To specify when and how often the scheduled task will be executed, click the calendar icon next to **Schedule**.

Change Test Schedule

Run at scheduled time(s)

Start at: 

☐ Randomize start times
Start at a random time within 30 minutes of the time specified above.

☒ Repeat task

Repeats:

Repeat every: week(s)

Repeat on: ☐ S ☒ M ☐ T ☐ W ☐ T ☐ F ☐ S

End:

Summary: At 2020-09-21 08:33:00 -07:00, repeat every 1 week(s) on Monday

Figure 8. Set or change test schedule

- a Next to **Start at** click the calendar icon and select the date and time to start the schedule.
- b Select **Randomize start times** to have the schedule start randomly within 30 minutes of the selected start time.
- c To set an interval for repeating the tests, select **Repeat task**.
- d Select **Hourly**, **Daily**, **Weekly**, or **Monthly** and set a repeat interval.
For **Weekly**, you can also select the days of the week for execution.
For **Monthly**, you can also select to repeat by day of the month or day of the week.

Using Test Services

Enable or disable a test schedule

- e Select an end date for the schedule. Select **Never** to continue the schedule without an end date. Select **On date** to specify a date to end the scheduled tasks.
 - f A summary of the selected options is displayed. Click **OK** to accept the selections.
- 6 When the schedule is complete, click **Save**.

Enable or disable a test schedule

To enable or disable a test schedule,

- 1 On the Test Services navigator, click the calendar icon to open the **Test Schedules** page.
- 2 In the test schedule list, select the test schedules to enable or disable.
- 3 At the top, click **Enable** to enable or **Disable** to disable the selected test schedules.

Change a test schedule

To change a test schedule,

- 1 Click the calendar icon to open the **Test Schedules** page.
- 2 From the list of test schedules, locate the test schedule you want to change, and click the **Edit** button on the right. 
- 3 Change the test schedule, as shown in **"Create a test schedule"** on page 29.

Delete a test schedule

To delete a test schedule,

- 1 Click the calendar icon to open the **Test Schedules** page.
- 2 From the list of test schedules, select the test schedules you want to delete, then at the top of the page, click **Delete**.
- 3 To confirm permanent deletion of the test schedule, click **Yes**.

Settings

Use the Settings tab to set up service user credentials and notifications.

Setting up credentials for running tests

For systems using authentication, configure the credentials for service users who will run tests. If no users are set in this section, any logged in user with correct privileges can start Test Services tests.

- 1 On the Test Services navigator, click the Settings icon. 
- 2 The Settings window opens.
- 3 Under Credential Settings, select and configure the user credentials for running Test Services tests:

Ad-hoc tests

Select the user credentials for running user-initiated tests.

Always use the following credentials for running user-initiated tests

Select this option to use the credentials you enter to run all user-initiated tests. If this box is not selected, the credentials of the currently logged in user are used for running user-initiated tests.

Scheduled tests

Select the credentials for running scheduled tests.

Always use the following credentials for running scheduled tests Select this option to use the credentials you enter to run scheduled tests. If this box is not selected, credentials of the user who created or modified the schedule are used for running scheduled tests.

- 4 Scroll to the bottom of the Settings page and click **Save**.

Setting up email notifications

You can set up Test Services to send a summary report email when verification tests are executed. You can configure notifications to:

- Send a daily summary email
- Send a weekly summary email
- Specify one or more recipients for email notifications

NOTE

Email is sent using OpenLab Shared Services. Make sure the email server is set up using the OpenLab Control Panel. If email is not configured in the OpenLab Control Panel, an error message appears, and you cannot set up notifications. For information on how to set up email in the OpenLab Control Panel, see the OpenLab Control Panel help.

- 1 On the Test Services navigator, click the Settings icon. 
- 2 The **Settings** window opens.
- 3 Scroll down to the Notification Settings. By default, notifications are disabled.
- 4 To enable notifications, under Notification Settings, click **Enable**. This makes the fields for setting up the notifications active.

Using Test Services

Setting up email notifications

Notification Settings
Configure global notification settings, applicable for current Test Services environment.

Notifications:
Enable or disable sending reports and notifications for test execution(s).

Send Emails To:
Enter one or more emails separated by ";". Based on the selections below, the system will send reports and notifications to these addresses.

Test Execution Report:

☐ Send Daily Test Execution Report at

☐ Send Weekly Test Execution Report at on

Summary of all completed tests and their results, executed during the previous 24 hours or 7 days from specified daily or weekly time. Specify the time to send the report in 'hh:mm' 24-hour format. For example, enter '6:30 PM' as '18:30'.

Immediate Notifications:

☐ Notify immediately on errors

☐ Notify immediately on warnings

Send the Test Execution Report immediately when test errors or warnings occur.

Notification Options:

☐ Exclude details for passed tests in the Test Execution Report

☐ Send no more than immediate notifications in minutes

Select if details for passed tests will be included in the Test Execution Report and limit the number of immediate notification emails sent.

Figure 9. Notification Settings

- 5 In the **Send Emails To** box, type one or more email addresses, separated by a semicolon. The system will send notifications to the email addresses specified, using the options selected in the following sections.
- 6 In the **Test Execution Report** section, select one or both:
 - **Send Daily Test Execution Report** - sends a daily report of all tests executed the previous day, including their results (ending at midnight)
 - **Send Weekly Test Execution Report** - sends a weekly summary report of all tests executed the previous week (ending on Sunday)
- 7 In the **Immediate Notifications** section, select one or both:
 - **Notify immediately on errors** - sends an email whenever an error is reported
 - **Notify immediately on warnings** - sends an email whenever a warning is reported
- 8 In the **Notification Options** section, select options for your notifications:

Using Test Services

Setting up email notifications

- **Exclude details for passed tests in the Execution Report** - select to include only details from failed tests
- **Send no more than** - select to limit the number of immediate notifications to a designated number of notifications within a specified number of minutes.

9 When you are finished, click **Save**.

Disabling notifications

To disable all email notifications,

- 1 On the Test Services navigator, click the Settings icon. 
The **Settings** window opens.
- 2 To disable notifications, under Notification Settings, click **Disable**. This makes the fields for setting up the notifications inactive.
- 3 Click **Save**.

Enabling notifications

To turn on notifications once they are disabled,

- 1 On the Test Services navigator, click the Settings icon. 
The **Settings** window opens.
- 2 To enable notifications, under Notification Settings, click **Enable**. This makes the fields for setting up the notifications, as well as the **Save** button, active. You can now set up new notifications or edit existing notifications.

Notification reports

Notification reports are emailed to the designated addresses. The reports include a summary of tests performed during the notification period, and include links to reports.

NOTE

If necessary, log in to the specified machine to view reports.

Using Test Services

Setting up email notifications

An example of a daily test execution report sent via notifications is shown in the following figure.

```
Test Services Daily Report from 3/30/2026 17:00 (+02:00) to 3/31/2026 17:00 (+02:00)

Server hostname: PERF28-WAD-SRV1

=== Summary by status ===

* Failed: 21
* Not Available: 12
* Aborted: 0
* Passed with additional info: 11
* Passed: 34

=== Failed (21) ===

* Machine: scl00051901a.germany.agilent.com (https://scl00051901a.germany.agilent.com:443/testservices)
* Test: OpenLab Software Installation Verification Test

* Machine: scl00051801a.germany.agilent.com (https://scl00051801a.germany.agilent.com:443/testservices)
* Test: OpenLab CDS Connectivity Test

* Machine: scl00051901c.germany.agilent.com
(https://scl00051901c.germany.agilent.com:52088/openlab/testservices)
* Test: OpenLab CDS Connectivity Test

* Machine: scl00051801c.germany.agilent.com
(https://scl00051801c.germany.agilent.com:52088/openlab/testservices)
* Test: OpenLab CDS Connectivity Test

* Machine: scl00051801c.germany.agilent.com
(https://scl00051801c.germany.agilent.com:52088/openlab/testservices)
* Test: OpenLab Software Installation Verification Test

* Machine: scl00051801a.germany.agilent.com (https://scl00051801a.germany.agilent.com:443/testservices)
* Test: OpenLab Software Installation Verification Test

* Machine: scl00051902a.germany.agilent.com (https://scl00051902a.germany.agilent.com:443/testservices)
* Test: OpenLab Software Installation Verification Test

* Machine: perf28-wad-srv1.germany.agilent.com (https://perf28-wad-srv1.germany.agilent.com:443/testservices)
```

Figure 10. Test Services daily report sent via notifications

In This Book

This document provides instructions on how to use Agilent Test Services.

www.agilent.com

© Agilent Technologies, Inc. 2022-2026

DocNo D0146323

04/2026

Published in USA

