



Agilent Lab Advisor

IT Administrator's Guide

Notices

Document Information

Document No: SD-29000258 Rev. F
Edition: 03/2024

Copyright

© Agilent Technologies, Inc.
2016-2024

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
Hewlett-Packard-Strasse 8
76337 Waldbronn, Germany

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.

Contents

1	Server Installation	4
	Server Installation	5
2	Network Licensing	6
3	Authentication	9
	Authentication Overview	10
	Setting Up Authentication	12
4	Agilent TCP Relay Service	15
	Agilent TCP Relay Service	16



1 Server Installation

This chapter describes how to install the Agilent Lab Advisor on a Windows Server 2016, 2019, or 2022 system.

Server Installation 5

Server Installation

You can install Lab Advisor on a Windows Server 2016, 2019, or 2022 system and make it available for use on client systems.

- 1 Install Lab Advisor on the Windows Server 2016, 2019, or 2022 PC as described in the section *Hard Drive Installation* of the *Agilent Lab Advisor User Manual*.
- 2 Install the role **Remote Desktop Services** on the Windows Server 2016, 2019, or 2022 server.
- 3 Publish the Lab Advisor installation on the Windows Server 2016, 2019, or 2022 server.
- 4 Create an RDP (Remote Desktop Protocol) file for the Lab Advisor installation.
- 5 Distribute the RDP file to all client systems.



2 Network Licensing

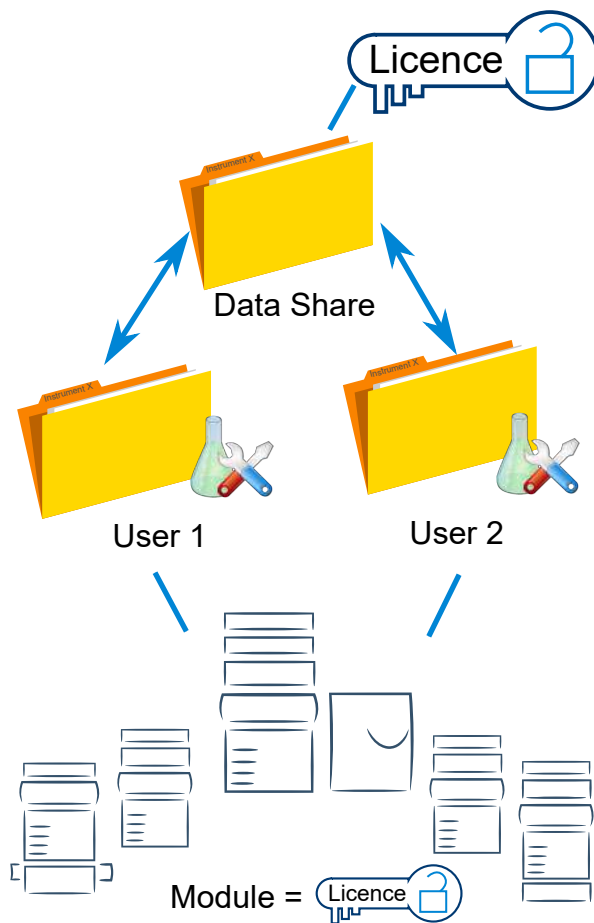
This chapter describes the licensing scheme for networked systems and gives instructions on how to set it up.

For Lab Advisor A.02.10 and later, the license scheme operates on the principle of one license per device, irrespective of the number of Lab Advisor installations. A single M8550A Advanced license in combination with add-on packs can be used simultaneously on multiple Lab Advisor installations. Licenses can be either typed in on any individual Lab Advisor installation or deployed via Data Sharing. In the case of Data Sharing, the licenses are stored in a network folder, and every Lab Advisor installation that has access to the folder can use the licenses.

With the 1-year advanced license M8556A (*per-user license*, licenses an unlimited number of hardware modules) Module License sharing is not appropriate and therefore not possible.

NOTE

Mixing *per-user* Unlimited licenses with the per module licenses will lead to unforeseen problems in conjunction with Data Sharing.



The basic requirements are one Advanced license (for the initial setup) with as many licenses as there are devices in the laboratory. All Lab Advisor installations with Basic licenses in the network will be automatically upgraded to the Advanced license when configuring the path to the Data Share..

Setting Up a Network License

- 1 In the network, create a folder that can be accessed by all Lab Advisor installations.
- 2 Install Lab Advisor on the first PC.
- 3 Start the Lab Advisor installation that you have just installed and navigate to **Configuration > Licenses**.
- 4 Type the *Advanced* License key into the **License Key** field and validate it by clicking **Add**.
Repeat the process for any Add-on license packs.
- 5 Navigate to **Configuration > Data Sharing**.
- 6 Click **Enable Data Sharing** and enter the URL of the network folder that you created in Step 1 into the Wizard.
- 7 In **Synchronize Folder**, select when you want to share instrument data in the network folder.
Select **Never** if you do not want to share instrument data.
- 8 Install Lab Advisor on all other PCs that you want to use,
- 9 On each PC, navigate to **Configuration > Data Sharing** and enable Data Sharing with the URL of the network folder that you created in Step 1 in **Data Share Folder**.
Network Licensing is now set up.



3 Authentication

This chapter describes the authentication options in Lab Advisor and gives instructions on how to set up authentication.

Authentication Overview 10

Lab Advisor Roles 11

Setting Up Authentication 12

Authentication Overview

The Advanced version of Lab Advisor provides two independent authentication providers:

- the **Legacy Lab Advisor Authentication Provider** is the equivalent of the **Traceability** authentication function of Lab advisor revisions prior to B.02.11.
- the **NT Authentication Provider** uses the authentication components of the Windows operating system.

When you enable authentication, you can select one or both providers to ensure that only authorized users have access to Lab Advisor functions and data.

NOTE

Before you enable Authentication, you must have set up and enabled at least one authentication provider with at least one user with the Administrator role, See the Lab Advisor online help for details.

Legacy Lab Advisor Authentication Provider

Users of the Legacy Lab Advisor Authentication Provider are set up during configuration using the **Add User** dialog box and are assigned one or more **Lab Advisor Roles** on page 11. During configuration, the administrator can specify that the user is required to log in with a password; this is equivalent to **Full Traceability** in Lab Advisor B.02.10 and earlier.

During configuration, the administrator decides whether to specify a password for each user or to allow users to specify their own passwords when they first log in to Lab Advisor.

All configured users and their details are shown in a table in the **Configuration** screen. The **Active** status of each user is denoted by a check box. This status is changed in the **Add User** dialog box.

NT Authentication Provider

Users of the NT Authentication Provider are those that are already configured as Windows users. During configuration, the administrator associates the **Roles** that are set up in Lab Advisor (see **Lab Advisor Roles** on page 11 with the **Groups** that are set up in the Windows networking environment. The administrator also decides whether or not users are allowed to edit their Lab Advisor profiles.

Lab Advisor Roles

A Lab Advisor role is characterized by a specified set of permissions. Lab Advisor roles are of two types:

- **Built-in** roles are provided by Agilent; the permissions can be viewed, but they cannot be edited or deleted. They include the **Administrator** role and the **Standard User** role.
- **User-defined** roles are set up by the administrator. User-defines roles have permissions that are specified by the administrator during configuration. They can be viewed and edited by users with Administrator permissions.

NOTE

The Administrator role must be assigned to at least one user in order to enable Authentication.

The **Roles** table in the **Authentication** screen shows the name of each role and its type.

Setting Up Authentication

Setting Up Data Sharing

If you intend to use Data Sharing, set it up *before* you set up Authentication.

1 Click **Configuration** in the **Tasks** list at the left of the Lab Advisor screen.

2 In the **Configuration** screen, click the **Data Sharing** icon.

The **Data Sharing** screen is displayed.

3 Click **Enable Data Sharing** and enter a URL to the folder that will contain the data from the Lab Advisor installation. Click **Next** and finish the wizard.

NOTE

Enabling or disabling a Data Share path has influence on the **Authentication** behavior. With Data Sharing, the **Authentication** configuration of the Data Share is synchronized. When Data Sharing is disabled the local **Authentication** configuration is used (when configured).

4 Restart Lab Advisor.

Data Sharing is activated and the uploaded data are available for review in the **Review Client**.

Setting Up Lab Advisor Roles

Lab Advisor provides two **Built-in** roles (**Administrator** and **Standard User**) with defined privileges. These roles cannot be edited or deleted. You can also set up additional **User-defined** roles with a different set of privileges.

1 Click **Configuration** in the **Tasks** list at the left of the Lab Advisor screen.

2 In the **Configuration** screen, click the **Authentication** icon.

The **Authentication** screen is displayed; the Roles are shown in a table in the lower half of the screen.

Select one of the roles from the table and click **View** to examine the privileges for that role.

3 To add a new role, click **Add**.

The dialog box is displayed to add a role.

Authentication

Setting Up Authentication

- 4 Enter a name for the new role in the **Role name** field.
- 5 Mark the check boxes for the privileges you want the new role to have.
- 6 When you are finished, click **Save**.
The new role is added to the **Roles** table, and is marked as **User-defined**.

Configuring the Provider

The two Authentication providers offered by Lab Advisor are configured differently.

- 1 In the **Authentication providers** section of the **Authentication** screen, select the provider you want to configure and click **Configure**.
The configuration dialog box for the selected provider is displayed.

Configuring the Legacy Lab Advisor Authentication Provider

- 1 To define users for the Legacy Lab Advisor Authentication Provider, click **Add** at the bottom right of the dialog box.
A dialog box is displayed to allow you to enter the user's details.
- 2 Provide the information for the user:
 - In the **General information** section, enter the details of the user.
 - Mark the **Allow user to set a new password on next login** to give the user the opportunity to change the Password specified in the **General Information** section at the next login.
 - Mark the **Active** check box to allow the user to log in. The state of this check box is shown in the **Configure Users** table
 - In the **Role membership** section, mark the check box(es) against the role(s) that you want the new user to belong to.
NOTE: You must assign at least one user to the **Administrator** role.
- 3 When you are finished, click **Save** to save the new user.
The new user is added to the **Configured Users** table. You can edit a user's details at any time by selecting the user in the table and clicking **Edit**.
- 4 In the **General Settings** section,

Authentication

Setting Up Authentication

- Mark the **Require users to enter a password to log in** check box to ensure that only authorized users with a Username and Password can use Lab Advisor. This is equivalent to **Full Traceability** in Lab Advisor revisions prior to B.02.11.
- Mark the **Allow users to change their own details and reset their password** check box to allow users to edit the entries in the Add User dialog box. If this check box is cleared, only users with administrator rights can make changes to users' details.

- 5 When configuration is complete, click the **X** button at the top right of the dialog box.

Configuring the NT Authentication Provider

- 1 In the **General Settings** section, mark the **Allow users to change their own details** check box to allow users to edit their entries in their Lab Advisor Profile when they are logged in to Lab Advisor.
- 2 In the **Group Settings** section, map Lab Advisor roles (built-in and user-defined) to the NT groups set up in the network. Type the name of the NT group in the appropriate field.
- 3 Click **Save settings** to validate the groups and save the associations.
Any NT groups that do not exist are marked in red as invalid; all invalid mappings must be corrected before you can save the mappings.
- 4 When configuration is complete, click the **X** button at the top right of the dialog box.

Enabling the Provider and enabling Authentication

- 1 In the **Authentication providers** section of the **Authentication** screen, select the provider(s) that you want to enable and click **Enable**.
- 2 Mark the **Enable authentication system** to enable the Lab Advisor authentication system.

NOTE

Before you mark this check box, ensure that you have set up and enabled at least one authentication provider with at least one user with the **Administrator** role,



4 Agilent TCP Relay Service

This chapter introduces the Agilent TCP Relay Service.

Agilent TCP Relay Service 16

Agilent TCP Relay Service

Agilent Lab Advisor revision B.02.09 and later is supplied with the Agilent TCP Relay Service, which replaces the Agilent Lab Advisor Relay Service. The TCP Relay Service has wider applicability than the Lab Advisor Relay Service, and provides a simplified user interface and password security. For full details of the Agilent TCP Relay Service, see the *Agilent TCP Relay Service Administrator's Guide*, available on the installation CD-ROM in the `\Support\RelayService` folder.

In This Book

This manual contains information for IT Administrators of the Agilent Lab Advisor.

The manual describes the following:

- Server Installation
- Network Licensing
- Authentication
- Agilent TCP Relay Service

www.agilent.com

© Agilent Technologies Inc. 2016-2024
Edition: 03/2024

Document No: SD-29000258 Rev. F

