

---

**README - Agilent Administration software**

---

**Contents**

**1. Introduction**

- 1.1 New Features with Agilent Administration software revision 1.0.1
- 1.2 New Features with Agilent Administration software revision 1.0.0

**2. System Configuration**

- 2.1 PC Hardware Requirements
- 2.2 Operating System Requirements

**3. Software Installation and Setup**

- 3.1 Installation Instructions
- 3.2 Uninstall Instructions

---

## 1 Introduction

---

This document provides last minute information and installation advice corresponding to the Agilent Administration software.

### 1.1 New Features with Agilent Administration Software revision 1.0.1

- Administration Software version 1.0.1 introduces compatibility with the ProteoAnalyzer system, to support compliance requirements.
- A white paper as resource for users of the Agilent ProteoAnalyzer system whose organizations must comply with US FDA Part 11 in Title 21 of the Code of Federal Regulations (CFR), and its EU analog, Eudralex Chapter 4, Annex 11 PN: 5994-8836EN is available.
- An updated version of the white paper as resource for users of the Agilent Fragment Analyzer system whose organizations must comply with US FDA Part 11 in Title 21 of the Code of Federal Regulations (CFR), and its EU analog, Eudralex Chapter 4, Annex 11 PN: 5994-8835EN is available.
- The Administration software allows users to set up accounts with a numerical username

### 1.2 New Features with Agilent Administration Software revision 1.0

- Includes workflow management, access control and user authentication.
  - Creation of projects, all analytical runs must occur within a project.
  - Assigning roles to users with the ability to create customized roles.
  - Creation of reports on projects, roles, users and system-wide activities.
  - Ability to set an electronic signature workflow.
- New elements are present within the Fragment Analyzer controller software and ProSize data analysis software. They are described in their respective User Manuals.
- A white paper as resource for users of the Agilent Fragment Analyzer system whose organizations must comply with US FDA Part 11 in Title 21 of the Code of Federal Regulations (CFR), and its EU analog, Eudralex Chapter 4, Annex 11, is being prepared.

---

## 2 System Configuration

---

### 2.1 PC Hardware Requirements

The Agilent Administration software tested requirements are as follows:

Processor: i5 or higher / **ARM processor not supported**

RAM: Physical: 8 GB

Display: Min resolution of 1280x1024 or 1280x800 pixels

Hard disk: 500 GB recommended

### 2.2 Operating System Requirements

The Agilent Administration software has been fully tested and is supported to run on

- Microsoft © Windows 10 64 bit with **English (US)** language settings
- and
- Microsoft © Windows 11 64 bit with **English (US)** language settings

The Agilent Administration software tested requirements are as follows:

Operating system:	Microsoft © Windows 10 64 bit 1803 or later
Operating system language:	English (US)
Operating system region settings:	English(US) in default configuration
Microsoft .NET:	Microsoft .NET:
Adobe Acrobat DC:	DC 19.012.0034 (English US) or higher required for reporting

OR

Operating system:	Microsoft © Windows 11 64 bit 24H2 (OS Build 26100.3775)
Operating system language:	English (US)
Operating system region settings:	English(US) in default configuration
Microsoft .NET:	4.6.2 or higher
Adobe Acrobat DC:	DC 19.012.0034 (English US) or higher required for reporting

---

## 3 Software Installation and Setup

---

### 3.1 Installation Instructions

The Administration software installation program 'Setup.exe' is a Microsoft Windows application. You must have Windows 11 including latest updates installed prior to installing the Administration software. To install the Administration software, perform the following steps:

1. Start Windows.
2. Before installing please make sure you have administrative rights to install the software on your computer. You need to have an account with **administrative rights** to successfully install all required components.
3. Close **all** running programs.
4. Download the installation program from the following link.
5. Select the Agilent Administration software 'Register for download' icon.
6. Complete the Personal Data information and select the 'Submit' icon.
7. Select the Agilent Administration software to download.
8. Open the 'Administration software.zip' to locate the 'Administration software.exe'.
9. Double left-click on the 'Administration software.exe'
10. If a Windows security message appears, select 'Yes', to allow the program to make changes on the computer.
11. When the window "The installer will install the Administration Software" appears, select "OK" to open a self-extracting zip file.
12. Select "I accept the license agreement" and "Next" when prompted to begin installation.
13. Re-start to ensure successful software installation.

### 3.2 Uninstall Instructions

If after installing the Administration software, you find it necessary to remove the software from your computer, perform the following steps:

Uninstall instructions for Windows 10:

1. From the settings panel, open 'System' and 'Apps & features'.
2. Select '**Agilent Administration software**' from the list of apps that can be uninstalled by Windows.
3. Click on the 'Uninstall' button to uninstall the Administration software.

Uninstall instructions for Windows 11:

1. From the settings panel, open 'Apps' and 'Installed apps'.
2. Select '**Agilent Administration software**' from the list of apps that can be uninstalled by Windows.
3. Click on the 'Uninstall' button to uninstall the Fragment Analyzer controller software.

If you encounter problems whilst upgrading the Administration software, please remove the existing Administration software installations. If this does not resolve the installation issues, please use the tool contained in the link below or set up the laptop with recovery media.

[http://support.microsoft.com/mats/Program\\_Install\\_and\\_Uninstall/](http://support.microsoft.com/mats/Program_Install_and_Uninstall/)

---

© Copyright 2025 Agilent Technologies, Inc. Use, reproduction, and distribution are subject to approval of Agilent Technologies.

---

End of Readme ...