

---

**README - Agilent Fragment Analyzer controller software**

---

**Contents**

**1. Introduction**

- 1.1 New Features with Fragment Analyzer controller software revision 5.0.0**
- 1.2 New Features with Fragment Analyzer controller software revision 4.0.0.11**
- 1.3 New Features with Fragment Analyzer controller software revision 4.0.0**
- 1.4 New Features with Fragment Analyzer controller software revision 3.1.0.12**

**2. System Configuration**

- 2.1 System Requirements**
- 2.2 PC Hardware Requirements**
- 2.3 Operating System Requirements**

**3. Software Installation and Setup**

- 3.2 Installation Instructions**
- 3.3 Uninstall Instructions**

---

## 1 Introduction

---

This document provides last minute information and installation advice corresponding to the Agilent Fragment Analyzer controller software.

### 1.1 New Features with Fragment Analyzer controller software revision 5.0.0

- Fragment Analyzer controller software to now work with new Administration software.
- Installer will be provided as two versions:
  - A combined package, consisting of Fragment Analyzer controller software, the new Administration software and the ProSize data analysis software, for users who need to meet compliance requirements (Fragment Analyzer Software Security Module).
  - A standalone Fragment Analyzer installer, for users who don't need to work under compliance regulations.
- Logins for Fragment Analyzer software Secure Module to be set using Administration software
- Simplified Hardware testing screen. Service-level functions no longer accessible on this screen.
- "Method Save Bug" fixed.
- Improved communication between instrument and PC.
- "Fill Capillaries with Dye" message has been removed.
- .PSDA file data extension produced as data acquisition from Fragment Analyzer controller software and readable with ProSize data analysis software when Fragment Analyzer Software Security Module is set up.
- Removed features:
  - Login features are removed for standalone version of the software.
  - Enhanced security features removed from controller software.
  - Removed the following menu options:
    - Configuration>Security Settings
    - Configuration>Email
    - Administration>User Maintenance
    - Administration>Change Password
    - Administration>Archive & Purge Database

### 1.2 New Features with Fragment Analyzer controller software revision 4.0.0.11

- Instrument chooser in configuration settings now displays the available Fragment Analyzer systems
- Windows 11 compatibility

### 1.3 New Features with Fragment Analyzer controller software revision 4.0.0

- Timestamps display in local time zone in Event and Error Log

### 1.4 New Features with Fragment Analyzer controller software revision 3.1.0.12

- Merged the 5400 Fragment Analyzer automated system into the 5200/5300 software
- Added solutions barcoding in the case that ZAG consumables want to be used. This is configurable.
- Added forced conditioning for ZAG consumables.
- Rebranded the software to Agilent name conventions. 5200 = 12,24 cap FA, 5300 = 48,96 cap FA, 5400 = Infinity
- Update to service password
- Added Agilent user
- EULA update and added to Help menu
- Agilent digital certificate added to final build
- Agilent serial number compatibility for device serial number

## 2 System configurations

---

### 2.2 System Requirements

#### 2.2.1 PC Hardware (minimum requirements)

The Fragment Analyzer controller software tested requirements are as follows:

Processor:	i5 or higher
RAM:	Physical: 8 GB
Display:	Min resolution of 1280x1024 or 1280x800 pixels
Hard disk:	500 GB recommended

#### 2.2.2 Operating System requirements

The Fragment Analyzer controller 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 Fragment Analyzer controller 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:	4.6.2 or higher
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

The Automatic reporting function in the Fragment Analyzer controller software requires Adobe Acrobat DC Reader version 19.012.0034 or later to be able to view and generate reports. All PCs supplied by Agilent Technologies will have this pre- installed. Reporting to .pdf format is the default.

---

### 3 Software Installation and Setup

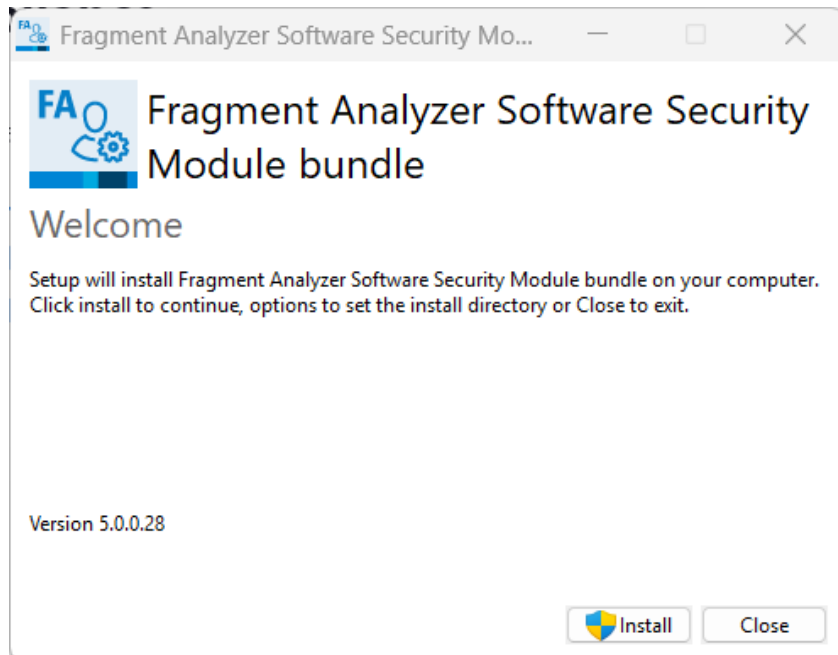
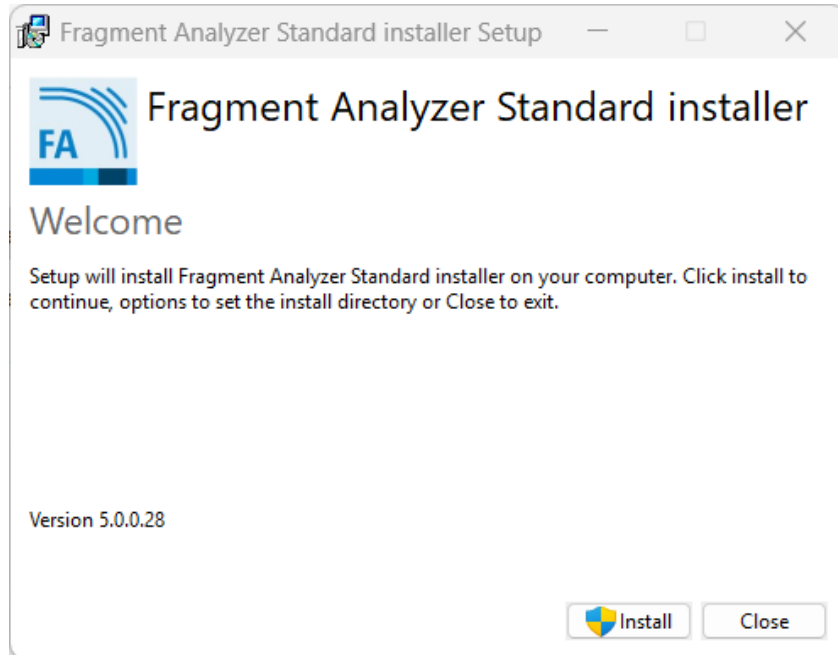
---

#### 3.2 Installation Instructions

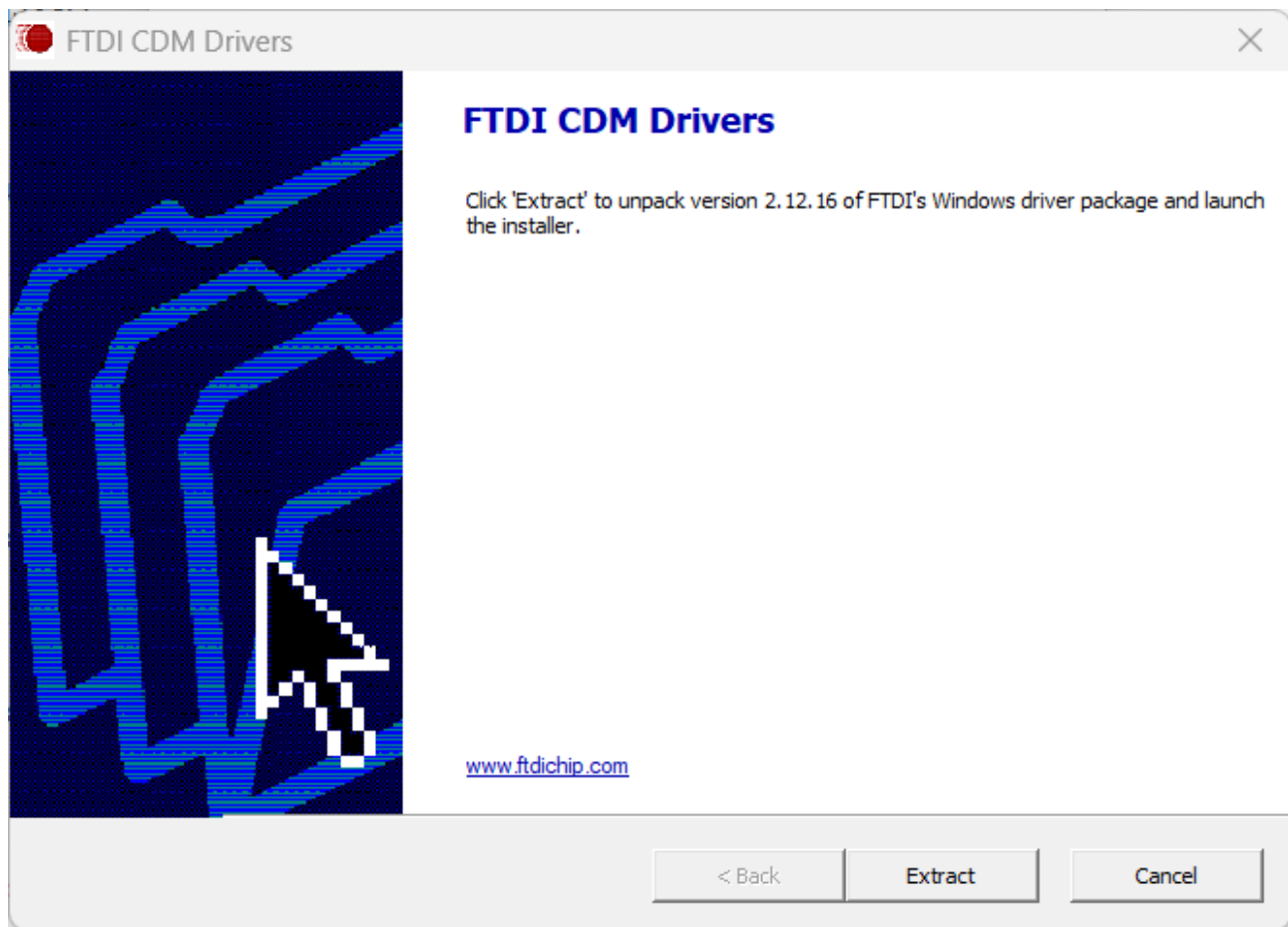
The Fragment Analyzer controller software installation program 'Fragment Analyzer Controller Software installer.exe' is a Microsoft Windows application. You must have Windows 10 or later installed prior to installing the Fragment Analyzer controller software.

To install the Fragment Analyzer controller 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, including all instances of any Fragment Analyzer controller software subprogram.
4. Download the installation program from the following link:  
<https://explore.agilent.com/Software-Download-Fragment-Analyzer-Prosize>
5. Select the proper installer to download. Either the Standard Installer or Fragment Analyzer Software Security Module bundle.
6. 'Download Revision 5.0.0' icon to download.
7. Open the 'Fragment Analyzer controller software installer and drivers V\_5.0.0' to locate the 'Fragment Analyzer controller software installer.exe'.
8. Double left-click on the 'Fragment Analyzer controller software installer.exe'.
9. If a Windows security message appears, select 'Yes', to allow the program to make changes on the computer.
10. When the following window appears select 'Install' to begin installation of the chosen software and drivers. Two starting screens shown below.

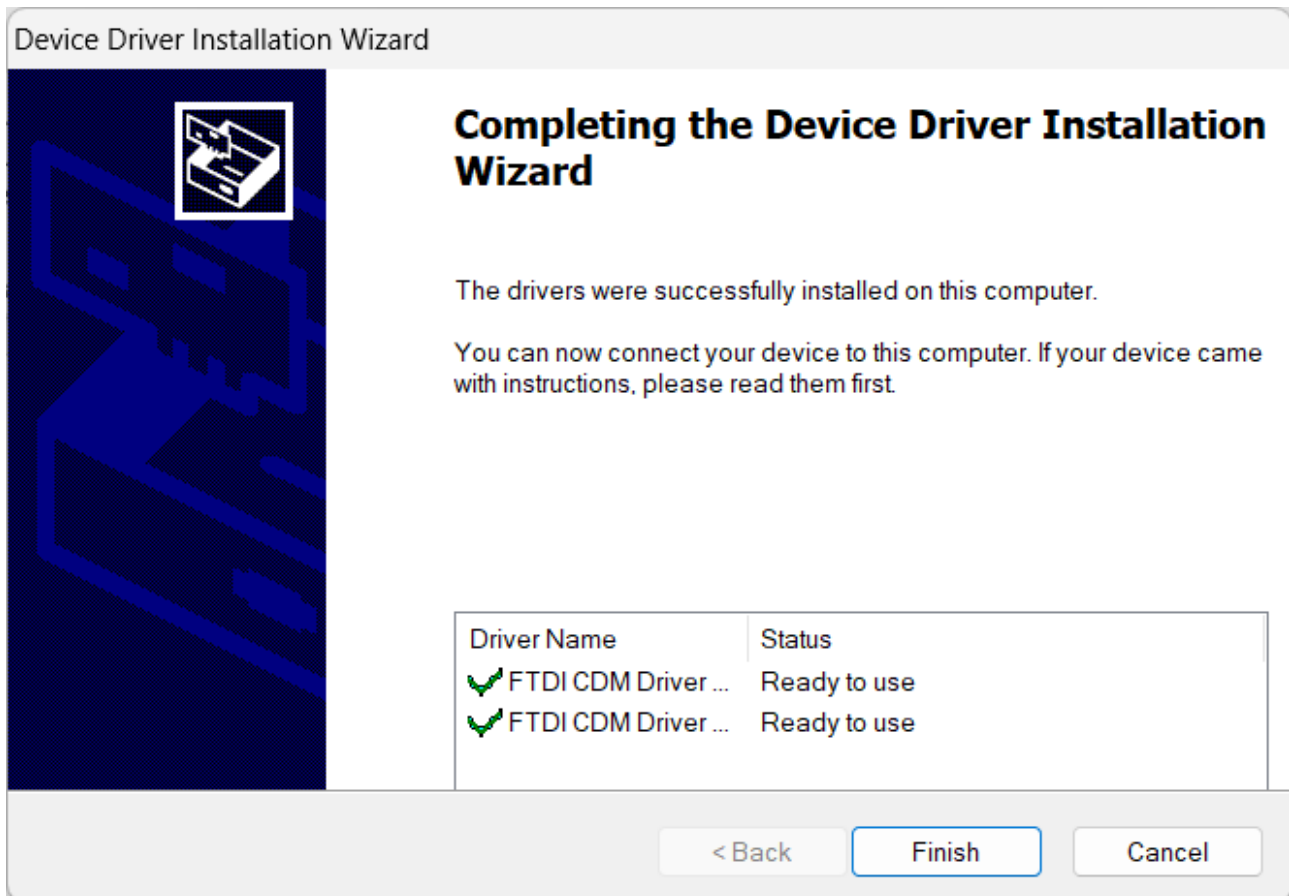


11. When the following window appears select 'Extract' to unpack the FTDI driver package.

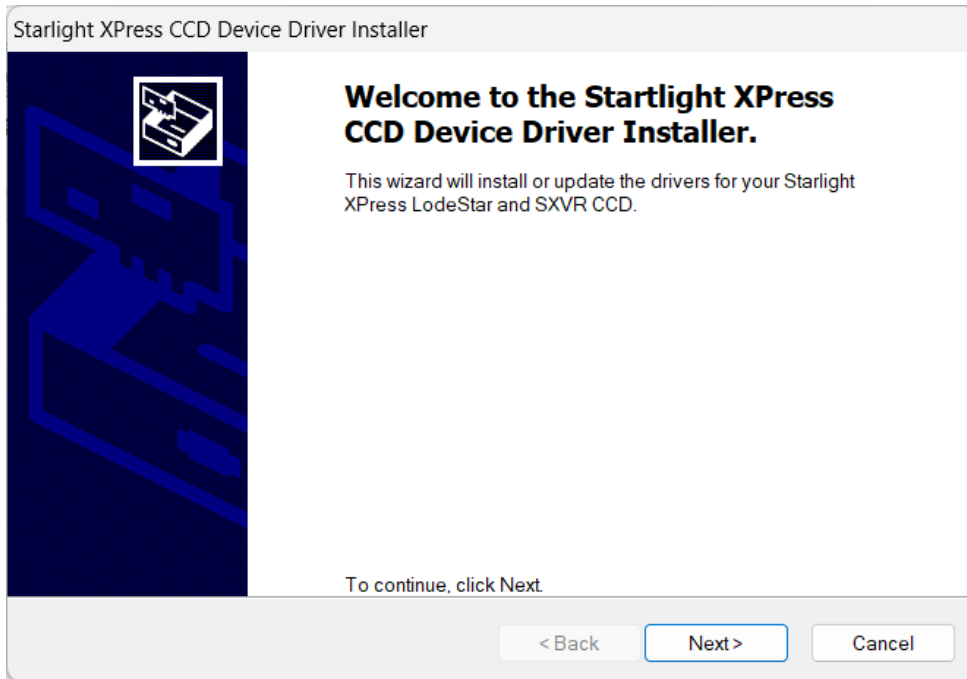


12. Select 'Next' to continue the installation, whenever needed.

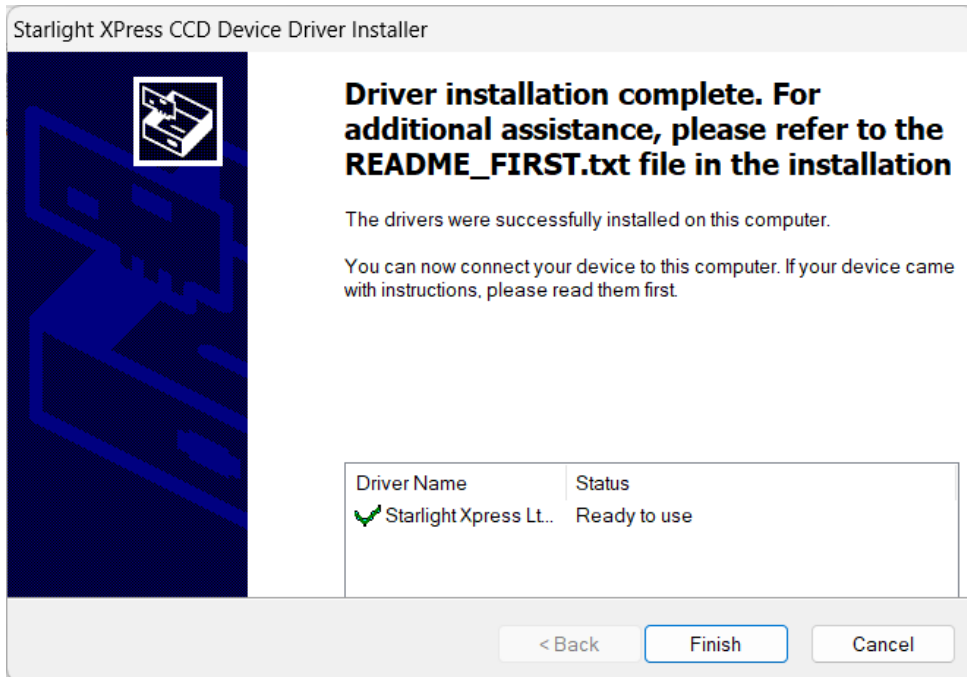
13. Once the FTDI drivers have been successfully installed, you will see the following window indicating the device driver is installed and is ready to use. Click 'Finish' to finish the installation.



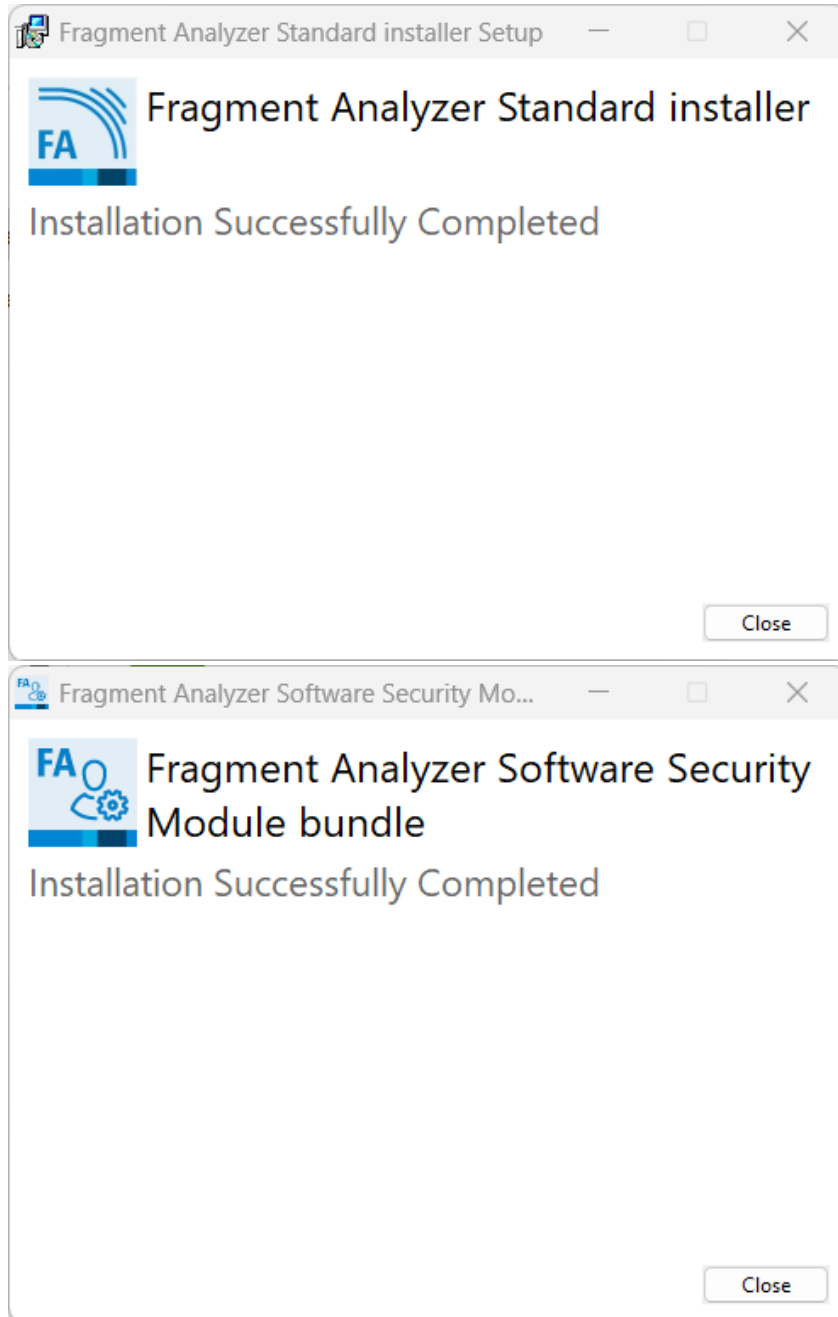
14. When the following window appears select 'Next' to begin installing the Starlight Xpress CCD Device Driver.



15. Once the Starlight Xpress CCD driver has been successfully installed, you will see the following window indicating the device driver is installed and is ready to use. Click 'Finish' to finish the installation.



16. Once the Fragment Analyzer controller software has been successfully installed, you will see the one of the following windows indicating the software is installed and is ready to use. Click 'Close' to finish the installation.



17. Re-start the PC to ensure successful software installation.

### 3.3 Uninstall Instructions

If after installing the Fragment Analyzer controller 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 '**Fragment Analyzer controller 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.

Uninstall instructions for Windows 11:

1. From the settings panel, open 'Apps' and 'Installed apps'.
2. Select '**Fragment Analyzer controller 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 Fragment Analyzer controller software, please remove the existing Fragment Analyzer controller 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/)

NOTE: If the Fragment Analyzer Software Security Module bundle was installed by accident, this security module must be uninstalled before attempting to run the Standard installer.

---

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

---

End of Readme ...