
README - Agilent ZAG DNA Analyzer controller software

Contents

1. Introduction
 - 1.1 New Features with ZAG DNA Analyzer controller software revision 2.2.0
 - 1.2 New Features with ZAG DNA Analyzer controller software revision 2.1.0.4
2. System Configuration
 - 2.1 System Requirements
 - 2.2 PC Hardware Requirements
 - 2.3 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 ZAG DNA Analyzer controller software.

1.1 New Features with ZAG DNA Analyzer controller software revision 2.2.0

- Improved communication between instrument and PC.

1.2 New Features with ZAG DNA Analyzer controller software revision 2.1.0.4

- Updated to newest libraries
- Rebranded to Agilent standards
- Updated service password.
- Added agilent user
- Removed ability to run a separation using perpetual barcodes.
- User input device serial number
- EULA update and added to Help menu
- Agilent digital certificate added to final build
- Windows 11 compatibility

2 System configurations

2.1 System Requirements

2.1.1 PC Hardware (minimum requirements)

The ZAG DNA Analyzer controller 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

2.1.2 Operating System requirements

The ZAG DNA 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 ZAG DNA 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

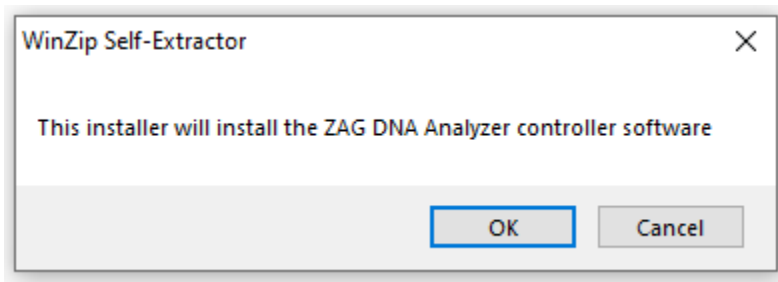
3 Software Installation and Setup

3.1 Installation Instructions

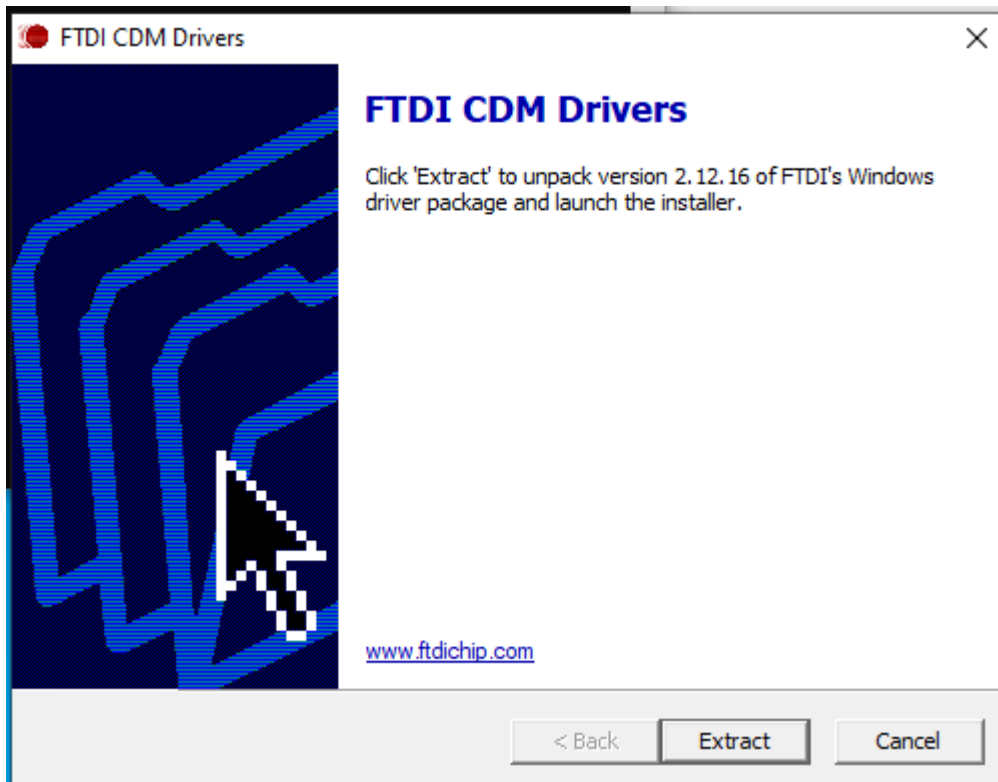
The ZAG DNA Analyzer controller software installation program 'ZAG DNA Analyzer controller software installer.exe' is a Microsoft Windows application. You must have Windows 10 or later installed prior to installing the ZAG DNA Analyzer controller software.

To install the ZAG DNA 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 ZAG DNA Analyzer controller software subprogram.
4. Download the installation program from the following link:
<https://explore.agilent.com/Software-Download-ZAG-DNA-Analyzer>
5. Select the 'Download Revision 2.2.0' icon to download.
6. Open the 'ZAG DNA Analyzer controller software installer and drivers V_2.2.0' to locate the 'ZAG DNA Analyzer controller software installer.exe'.
7. Double left-click on the 'ZAG DNA Analyzer controller software installer.exe'.
8. If a Windows security message appears, select 'Yes', to allow the program to make changes on the computer.
9. When the following window appears select 'OK' to open a self extracting zip file.

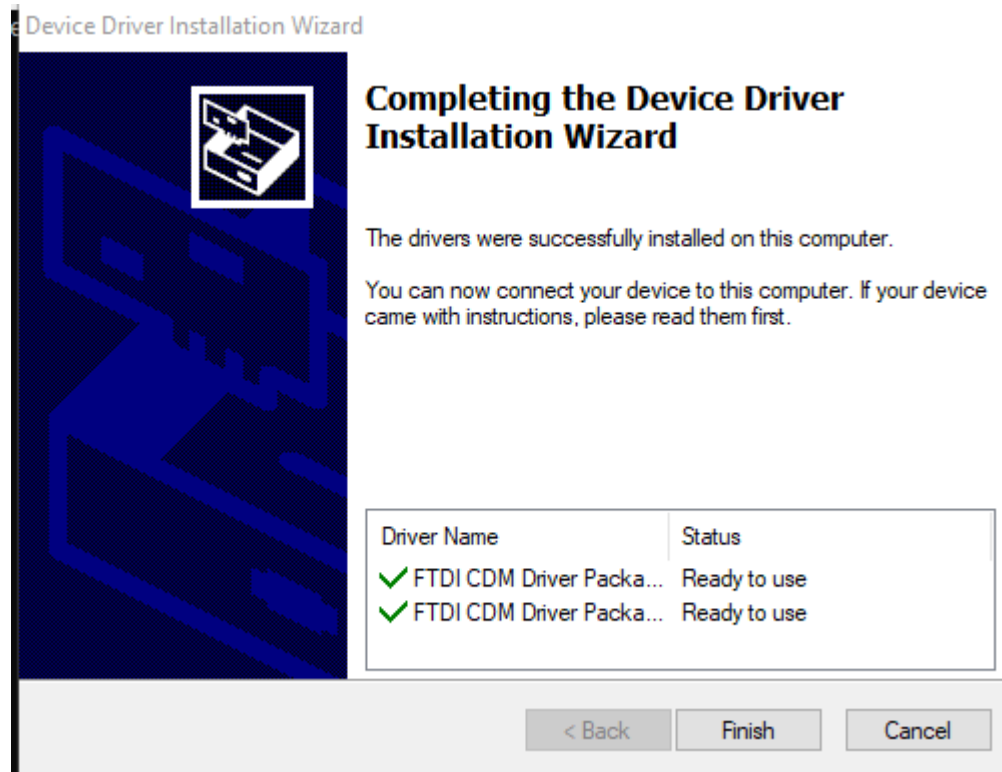


10. When the following window appears select 'Extract' to unpack the FTDI driver package.

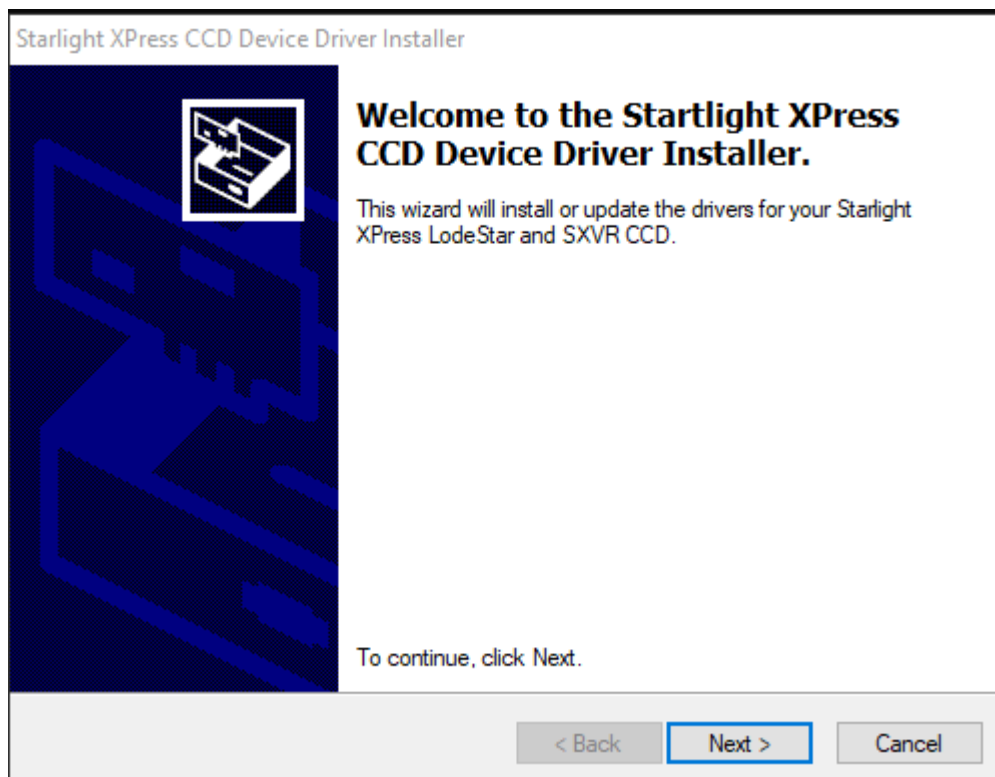


11. Select 'Next' to continue the installation, whenever needed.

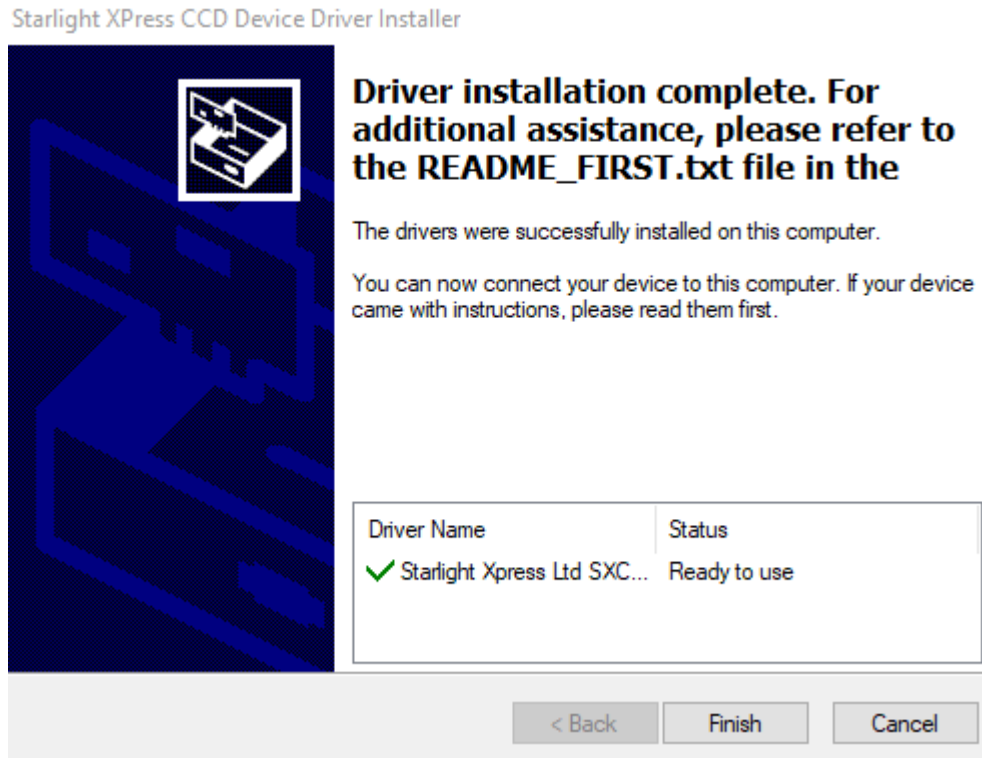
12. Once the FTDI drivers have been successfully installed, you will see the following window indicating the device driver is installed and is ready to use. Click 'Finish' to finish the installation.



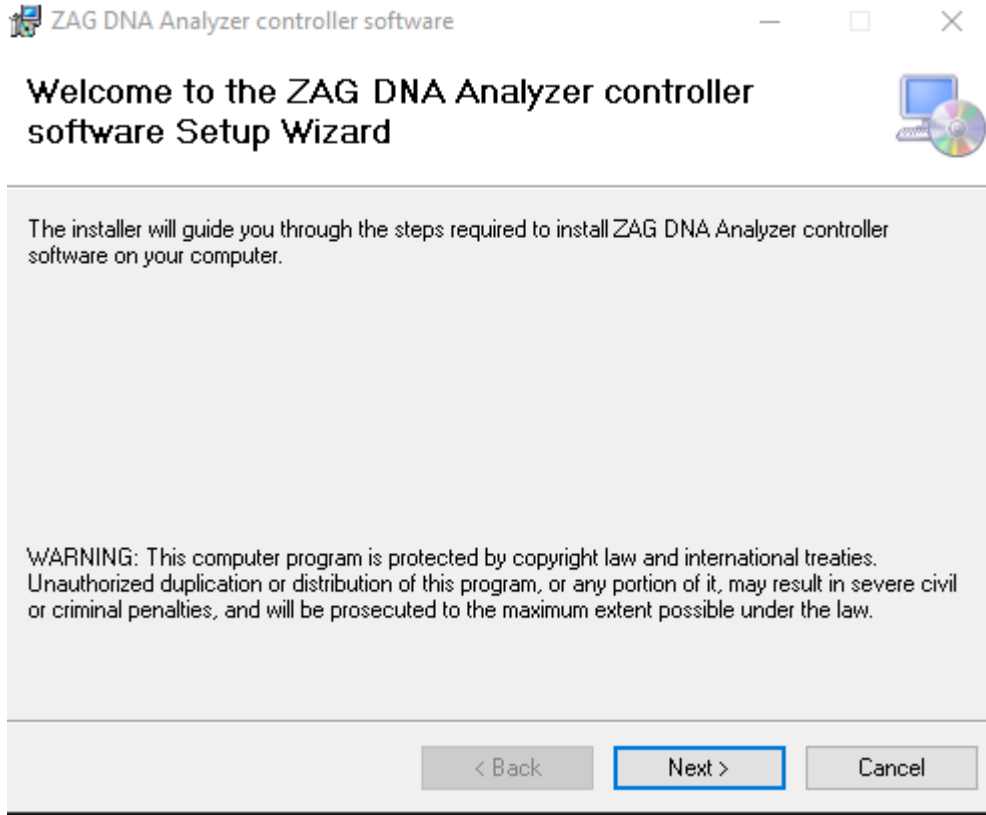
13. When the following window appears select 'Next' to begin installing the Starlight Xpress CCD Device Driver.



14. 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.

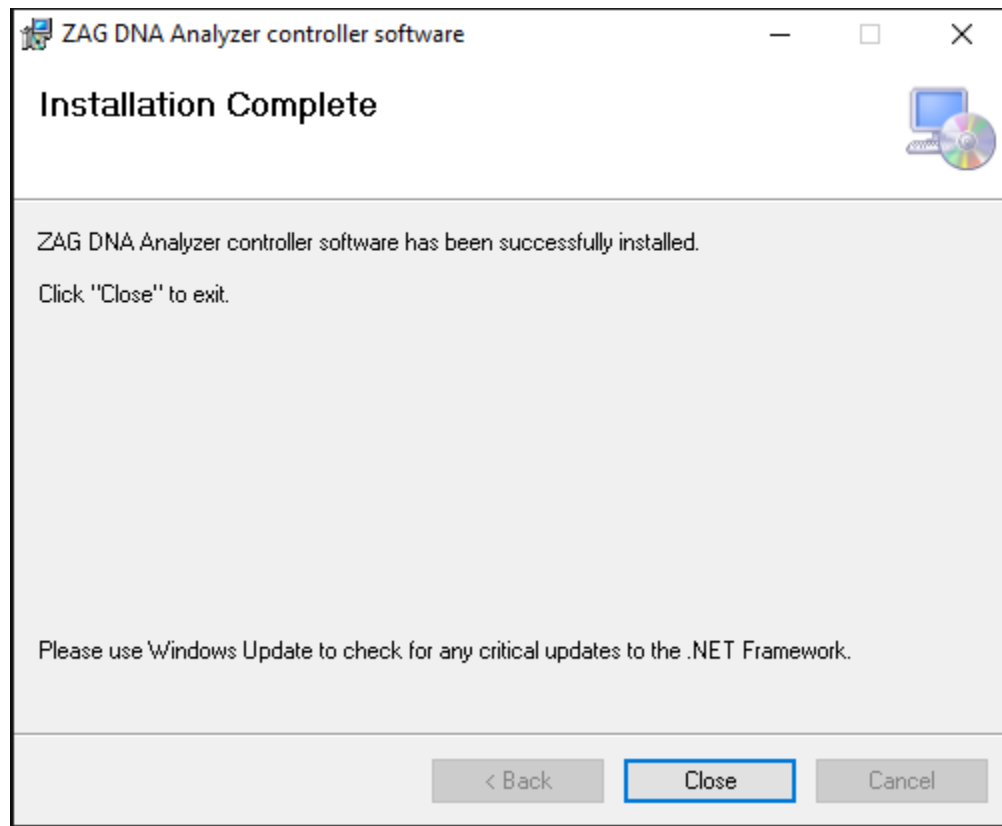


15. When the following window appears select 'Next' to begin installing the ZAG DNA Analyzer controller software.



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

17. Once the ZAG DNA Analyzer controller software has been successfully installed, you will see the following window indicating the software is installed and is ready to use. Click 'Close' to finish the installation.



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

3.2 Uninstall Instructions

If after installing the ZAG DNA 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 '**ZAG DNA Analyzer controller software**' from the list of apps that can be uninstalled by Windows.
3. Click on the 'Uninstall' button to uninstall the ZAG DNA Analyzer controller software.

Uninstall instructions for Windows 11:

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

If you encounter problems whilst upgrading the ZAG DNA Analyzer controller software, please remove the existing ZAG DNA 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/

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

End of Readme ...