
README - Agilent Femto Pulse controller software

Contents

1. Introduction

1.1 New Features with Femto Pulse controller software revision 2.0.0.3

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 Femto Pulse controller software.

1.1 New Features with Femto Pulse controller software revision 2.0.0.3

- Updated to Agilent branding standards
- Updated service password
- Added agilent user
- Agilent serial number compatibility for device serial number
- EULA update and added to Help menu
- Agilent digital certificate added to final build
- Windows 11 compatibility

2 System configurations

2.2 System Requirements

2.2.1 PC Hardware (minimum requirements)

The Femto Pulse controller software tested requirements are as follows:

Processor:	i5 or higher
RAM:	Physical: Minimum 4 GB for Windows 10 Minimum 8 GB for Windows 11
Display:	Min resolution of 1280x1024 or 1280x800 pixels
Hard disk:	500 GB recommended

2.2.2 Operating System requirements

The Femto Pulse 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 Femto Pulse 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 Femto Pulse 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.

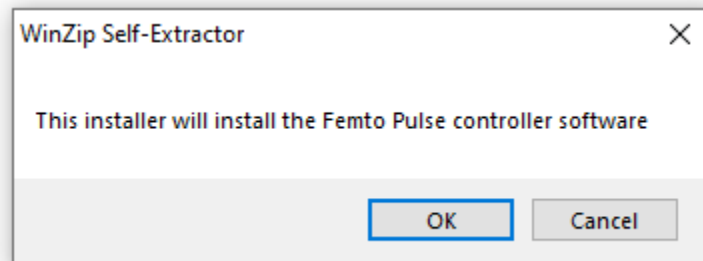
3 Software Installation and Setup

3.1 Installation Instructions

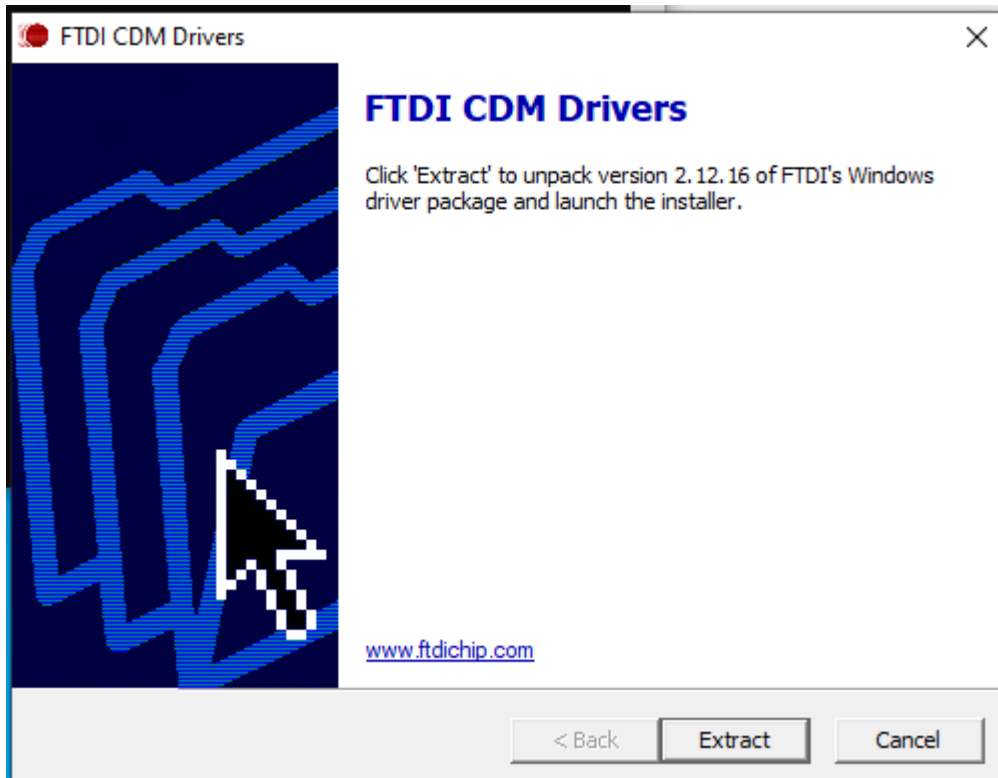
The Femto Pulse controller software installation program 'Femto Pulse Controller Software installer.exe' is a Microsoft Windows application. You must have Windows 10+ including latest updates installed prior to installing the Femto Pulse controller software.

To install the Femto Pulse 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 Femto Pulse controller software subprogram.
4. Download the installation program from the following link:
<https://explore.agilent.com/Software-Download-Femto-Pulse-Prosize>
5. Select the 'Download Revision 2.0.0.3 (35MB)' icon to download.
6. Open the 'Femto Pulse controller software installer and drivers V_2.0.0.3' to locate the 'Femto Pulse controller software installer.exe'.
7. Double left-click on the 'Femto Pulse 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.

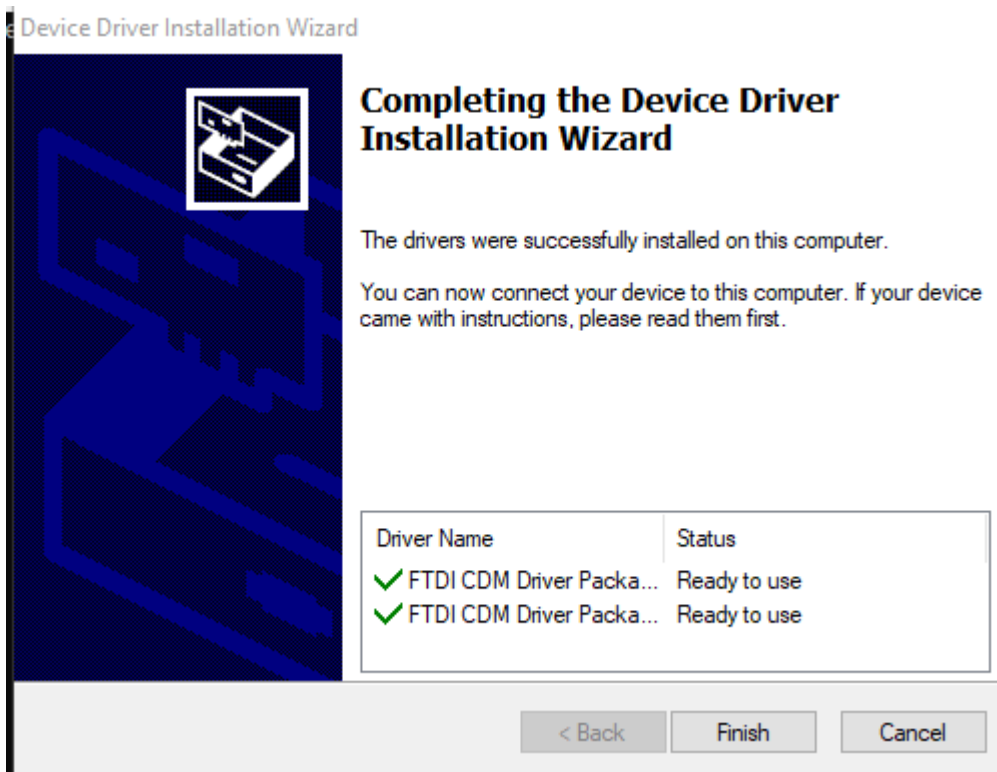


10. When the following window appears select 'Extract' to unpack the FTDI driver package.

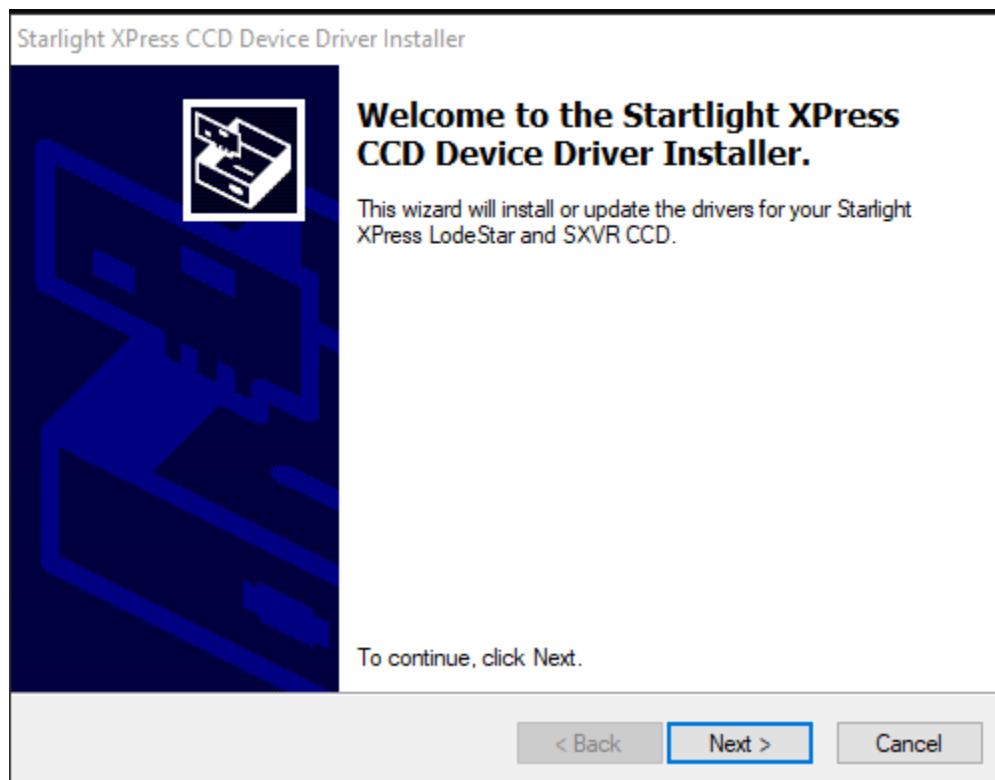


11. Select 'Next' to continue the installation, whenever needed.

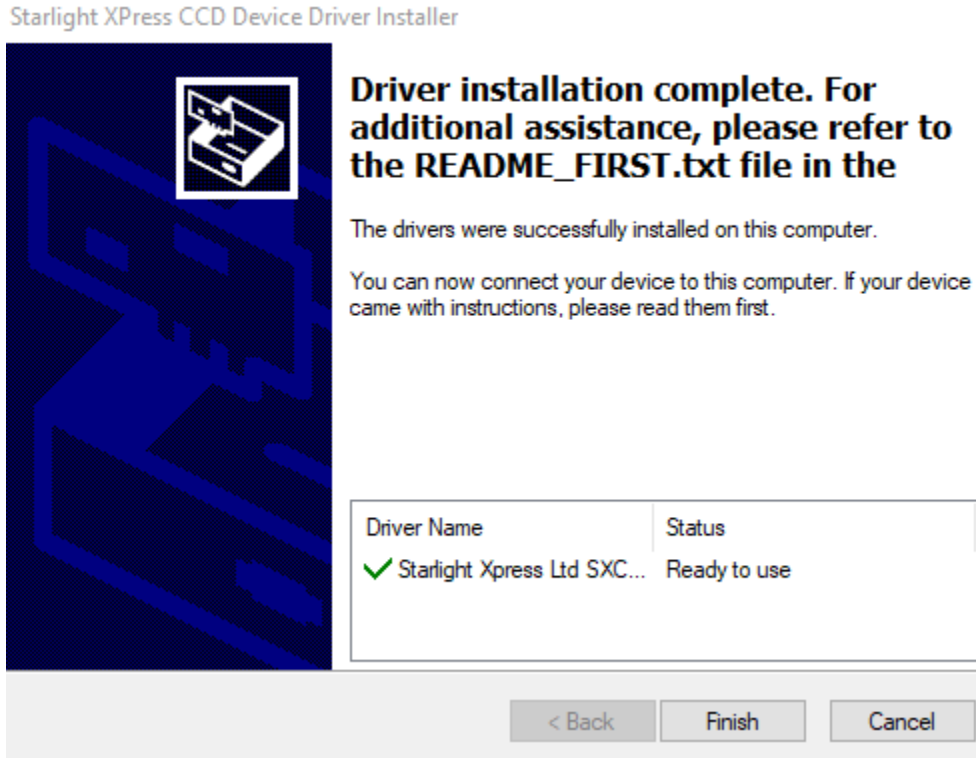
12. Once the FTDI drivers have been successfully installed, you will see the following window indicating the device driver is installed and is ready to use. Click 'Finish' to finish the installation.



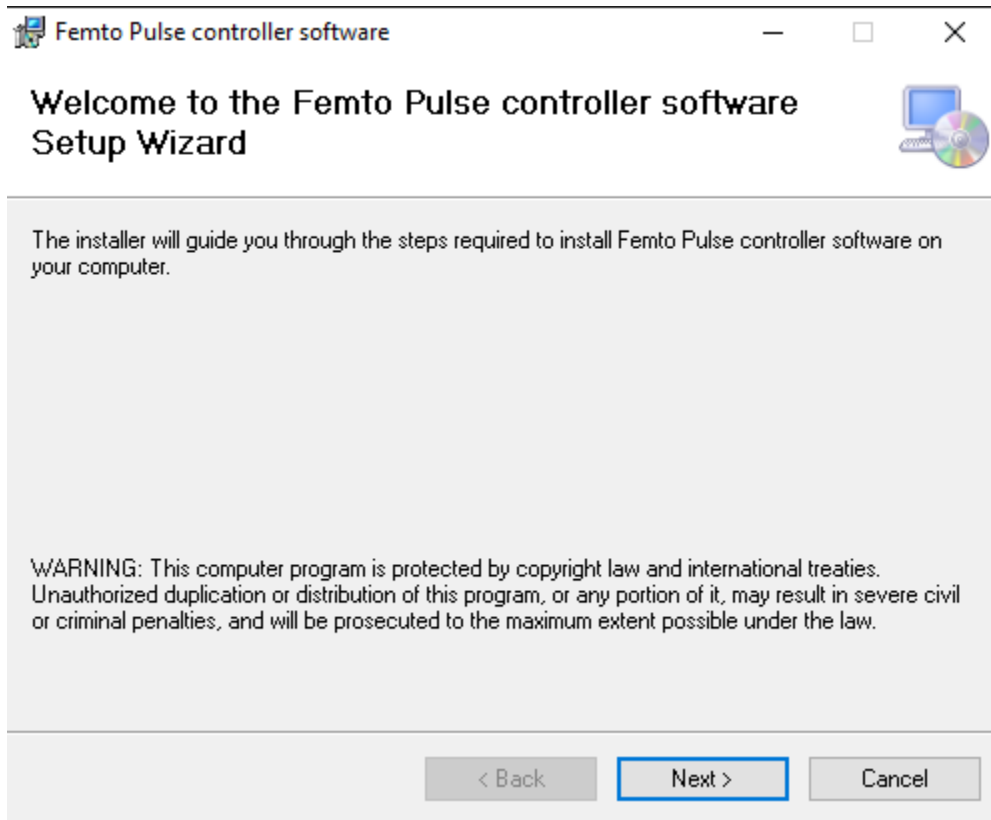
13. When the following window appears select 'Next' to begin installing the Starlight Xpress CCD Device Driver.



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.

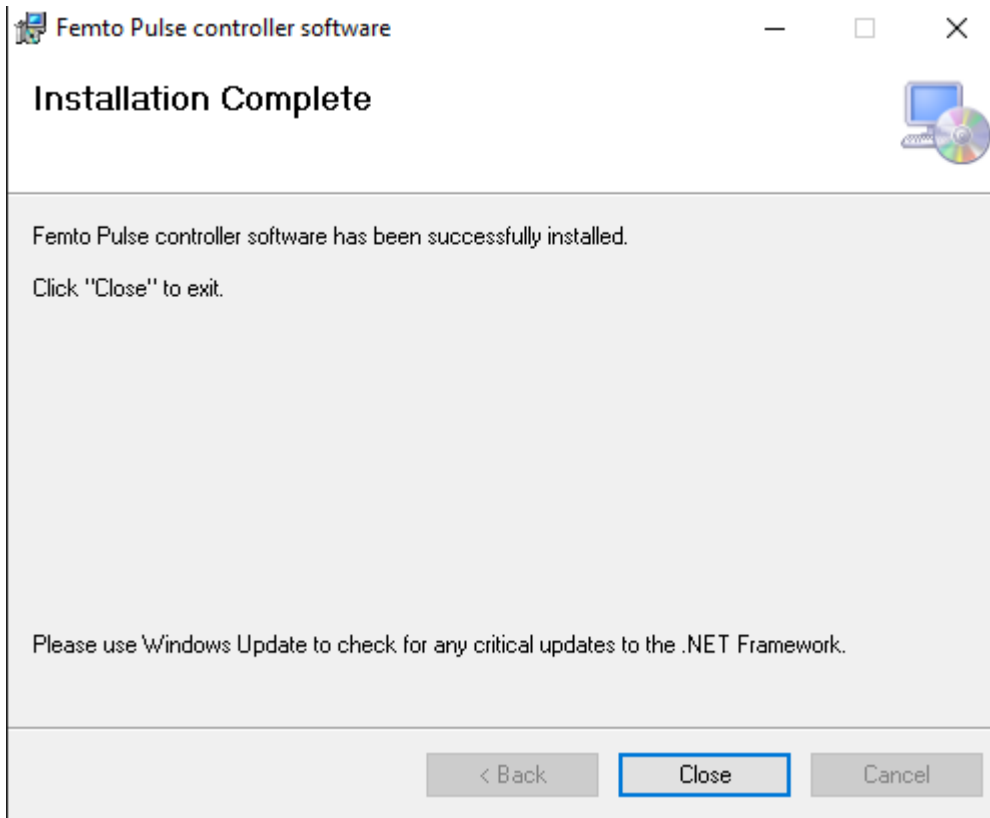


15. When the following window appears select 'Next' to begin installing the Femto Pulse controller software.



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

17. Once the Femto Pulse 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 Femto Pulse 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 '**Femto Pulse controller software**' from the list of apps that can be uninstalled by Windows.
3. Click on the 'Uninstall' button to uninstall the Femto Pulse controller software.

Uninstall instructions for Windows 11:

1. From the settings panel, open 'Apps' and 'Installed apps'.
2. Select '**Femto Pulse controller software**' from the list of apps that can be uninstalled by Windows.

3. Click on the 'Uninstall' button to uninstall the Femto Pulse controller software.

If you encounter problems whilst upgrading the Femto Pulse controller software, please remove the existing Femto Pulse 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 2021 Agilent Technologies, Inc. Use, reproduction, and distribution are subject to approval of Agilent Technologies.

End of Readme ...