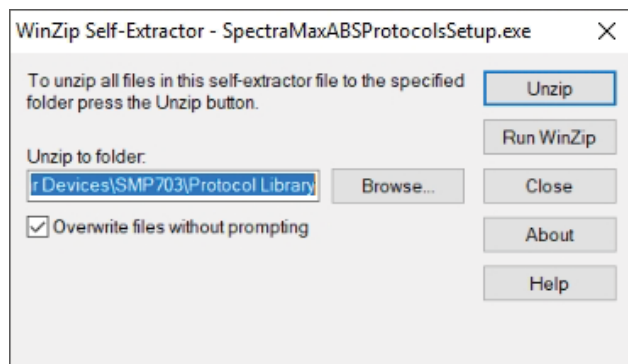


- In the WinZip Self-Extractor - SpectraMaxABSProtocolsSetup.exe dialog confirm that the **Unzip To Folder** field displays **C:\ProgramData\Molecular Devices\SMP703\Protocol Library**. If not, click **Browse** and navigate to and select the **C:\ProgramData\Molecular Devices\SMP703\Protocol Library** folder.



- In the WinZip Self-Extractor confirmation message, click **OK**.
- In the WinZip Self-Extractor - SpectraMaxABSProtocolsSetup.exe dialog, click **Close**.
- Before you start the software, power on the instrument you connect to the computer.

Get Help

For more information about how to use the SoftMax Pro Software, see the SoftMax Pro Software application help or user guide. To view the user guides, use a PDF viewer such as Adobe Reader. The software user guide, the formula reference guide, and user guides for the supported instruments can be found in the Windows Start menu:

Start > All Programs > Molecular Devices > SoftMax Pro 7.0.3

SpectraMax® ABS Reader Plug-in Installation Guide for SoftMax® Pro Software 7.0.3

Thank you for purchasing the SpectraMax ABS or SpectraMax® ABS Plus Microplate Reader. You operate this instrument from a computer that runs the SoftMax Pro Data Acquisition and Analysis Software version 7.0.3 with the SpectraMax ABS Reader Plug-in package installed. This document provides instructions for how to install the SpectraMax ABS Reader Plug-in package after you install the SoftMax Pro Software version 7.0.3 on the computer.

You should install the SoftMax Pro Software on the computer before you set up the instrument. Please be aware that some updates to the SoftMax Pro Software require a purchase. Contact Molecular Devices before you update your software. To download the SoftMax Pro Software version 7.0.3 and the SpectraMax ABS Reader Plug-in package, visit: www.MolecularDevices.com/SMPdownloadsite.

The SpectraMax ABS Reader Plug-in package updates the SoftMax Pro Software to support the operation of the SpectraMax ABS and SpectraMax ABS Plus Microplate Readers along with related online help instructions. The update also includes the default protocols and the validation protocols for the SpectraMax ABS and SpectraMax ABS Plus Microplate Readers.

You must install SoftMax Pro Software version 7.0.3 on the computer before you can install the SpectraMax ABS Reader Plug-in package. See the *SoftMax Pro Data Acquisition and Analysis Software Installation Guide* for computer system requirements and installation instructions.

Install the SpectraMax ABS Reader Plug-in

Before you can install the SpectraMax ABS Reader Plug-in you must install the SoftMax Pro Software version 7.0.3.

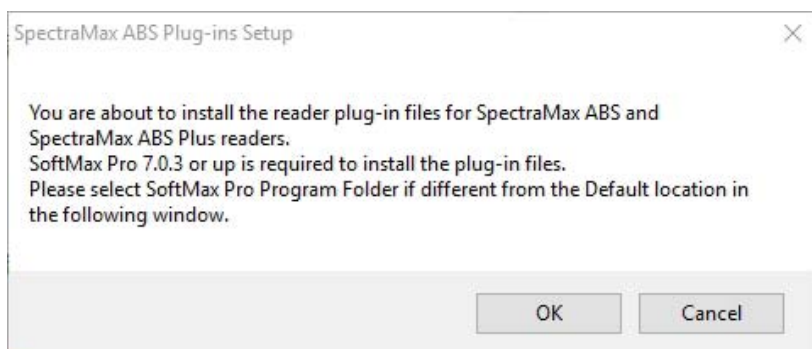
If you have not yet installed the SoftMax Pro Software on your computer, follow the installation instructions included with the software installation package.



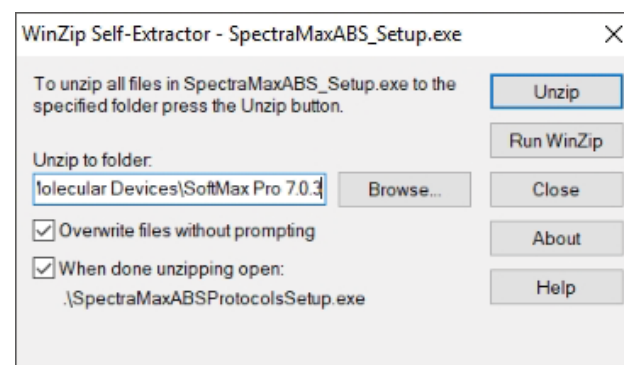
Note: You should disable the anti-virus program on the computer before you install SoftMax Pro Software, as it might interfere with the installation process.

To install the SpectraMax ABS Reader Plug-in:

1. Power off all connected instruments and close all versions of the SoftMax Pro Software on the computer.
2. Download or copy the SpectraMax ABS Reader Plug-in package from www.MolecularDevices.com/SMPdownloadsite to the computer running the SoftMax Pro Software.
3. Double-click the **SpectraMaxABS_Setup.exe** file to start the installation program.
4. When the SpectraMax Plug-ins Setup dialog displays, read the message and click **OK**.



5. In the WinZip Self-Extractor - SpectraMaxABS_Setup.exe dialog, confirm that the **Unzip To Folder** field displays C:\Program Files (x86)\Molecular Devices\SoftMax Pro 7.0.3. If not, click **Browse** and navigate to and select the **C:\Program Files (x86)\Molecular Devices\SoftMax Pro 7.0.3** folder.



6. Leave the two check boxes selected and click **Unzip**.
7. A message displays to indicate three files were successfully unzipped. Click **OK**.
8. In the SpectraMax ABS Protocol Setup dialog, click **OK**.

