

# Getting Started (Installing Protocols)

The *SpectraTest® ABS2 Absorbance Validation Plate User Guide* and the protocols for the SpectraTest ABS2 Absorbance Validation Plate are on the USB drive that shipped with your validation plate. After you install the protocols, follow the instructions in the *SpectraTest ABS2 Absorbance Validation Plate User Guide* to enter certificate information into the protocols and perform the validation tests.

This document provides instructions for how to install the protocols into the SoftMax® Pro Data Acquisition and Analysis Software Protocol Library.

## Installing the Protocols

The SpectraTest ABS2 Absorbance Validation Plate protocols included on the USB drive are compatible with the SoftMax Pro Software versions 6.5.1, 7.0.3, 7.1, and higher.



**Note:** SoftMax Pro Software version 7.1.1 is for the GxP edition only.

If needed, you can obtain the latest version of the validation protocols by contacting Molecular Devices support via the web site <https://www.moleculardevices.com/support.html>

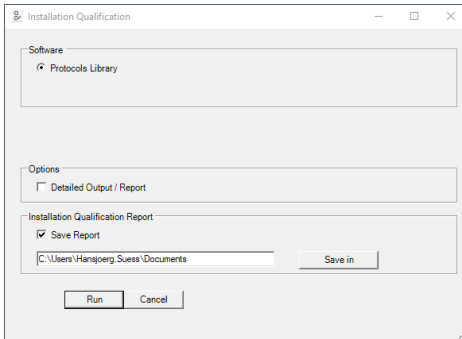
To install the SpectraTest ABS2 Absorbance Validation Plate protocols:

1. Power off all connected instruments and close all versions of the SoftMax Pro Software on the computer.
2. Insert the USB drive that shipped with the validation plate into a USB slot in the computer and open Windows Explorer.
3. Locate the .zip file that correspond to the version of the SoftMax Pro Software you are running and unzip it to extract the files.
4. Locate the executable file that corresponds to the version of the SoftMax Pro Software you are running and double-click it to start the installation program.
5. When the "Do you want to allow this app to make changes to your device?" message displays, click **Yes**.
6. When the System Updater dialog displays, you can click **See Log** for additional information and then click **Install**.



## SpectraTest ABS2 Protocol Installation

- In the Installation Qualification dialog, you can run a report that includes what was installed for GxP purposes.

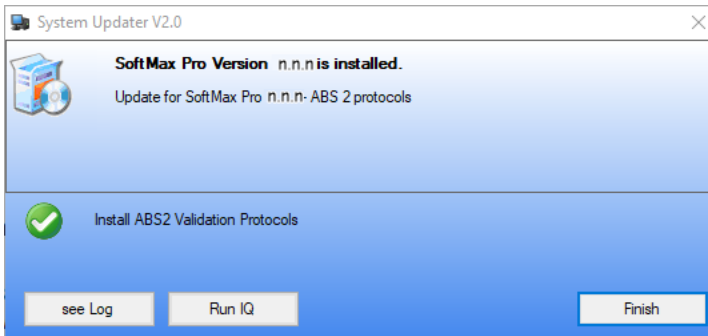


- Select the **Detailed Output/Report** check box to create an installation report that includes a list of the files that were added to the protocol library. Clear the **Detailed Output/Report** check box to create an installation report with a confirmation that files were added to the Protocol Library.
- Select the **Save Report** check box to define the location to save the report and to save the report for future reference. Clear the **Save Report** check box to not save the report for future reference.
- Click **Run** to run the report or click **Cancel** if you do not want to run the report.



**Note:** You can restart the System Updater at any time and click **Run IQ** to run the report in the future.


- After the installation completes, you can click **See Log** for a status message
- In the System Updater dialog, click **Finish**.

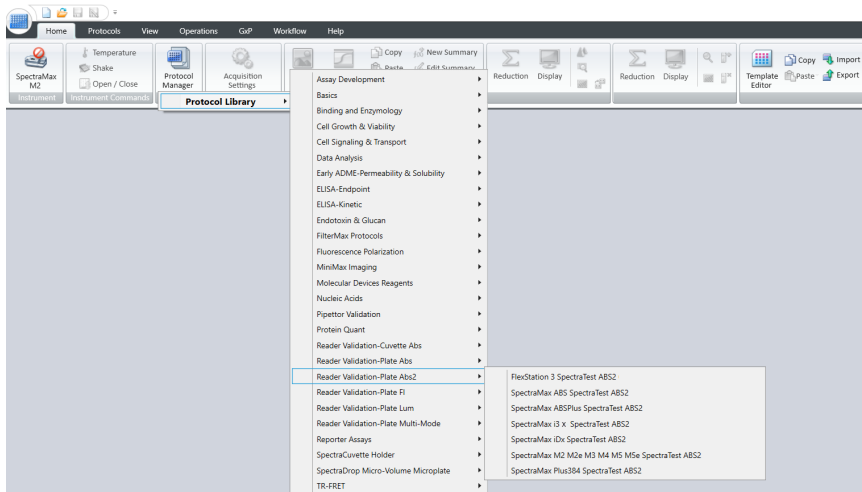


- Leave the USB drive in the computer and leave Windows Explorer open.
  - The instructions for how to use the SpectraTest ABS2 Absorbance Validation Plate are contained in the SpectraTest\_ABS2\_UserGuide.pdf file.
  - The information required to certify the protocols is contained in the (SN\_Date\_Cert#)\_EZinCert.pdf file.

## Verifying Installation of the Protocols

To verify installation of the protocols:

1. Start the SoftMax Pro Software.
2. Click the  **Protocol Manager** icon, expand the **Protocol Library**, and note that a subfolder named **Reader Validation-Plate Abs2** exists in the Protocol Library. The protocols in the Reader Validation-Plate Abs2 folder are dependent on the version of the SoftMax Pro Software you are running.



The SpectraTest ABS2 protocols are compatible with SoftMax Pro Software versions 7.1 and higher. SoftMax Pro Software version 7.1.1 is GxP edition only.

The following table lists protocol compatibility with earlier software versions.

**Table 1-1: Protocols**

Protocol Name	Version 6.5.1	Version 7.0.3
FlexStation 3 SpectraTest ABS2		✓
SpectraMax ABS SpectraTest ABS2		✓*
SpectraMax ABSPlus SpectraTest ABS2		✓*
SpectraMax i3(x) SpectraTest ABS2	✓	✓
SpectraMax iDx SpectraTest ABS2		✓
SpectraMax M2 M2e M3 M4 M5 M5e SpectraTest ABS2	✓	✓
SpectraMax Plus384 SpectraTest ABS2	✓	✓

\* Requires the installation of the SpectraMax ABS Reader plug-in that you probably installed if you are using a SpectraMax ABS microplate spectrophotometer.

## Get Help

For more information about how to use the SoftMax Pro Software, see the *SoftMax Pro Data Acquisition and Analysis Software User Guide* or the application help.

The SoftMax Pro Software installation places a copy of the microplate reader user guides (.pdf files) in the following location on the computer:

C:\ProgramData\Molecular Devices\User Guides

The most up-to-date microplate reader user guides are available on the Molecular Devices Knowledge Base at:

[support.moleculardevices.com](http://support.moleculardevices.com).

To access the Knowledge Base from the software:

- Select the **Home** tab, click **Contact Us**, and then select **Knowledge Base**.

### **For research use only. Not for use in diagnostic procedures.**

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/patents>

Product manufactured by Molecular Devices, LLC.  
3860 N. First Street, San Jose, California, 95134, United States of America.  
Molecular Devices, LLC is ISO 9001 registered.  
©2022 Molecular Devices, LLC.  
All rights reserved.