

Installing and configuring Sutter Instruments Lambda filter wheel\shutter devices

ABSTRACT

These instructions are for the installation of the Sutter instruments Lambda units under Windows 2000 professional and Windows XP professional only.

NOTES

This procedure applies to Meta Imaging Series software versions 6.0rx to 6.3rx.

OVERVIEW

HARDWARE SETUP AND TEST

Complete the following procedure to setup the filter wheel and/or shutter for use with your Meta Imaging Series software.

1. Turn the power switch off and plug the power cord in to the power connector located at the rear of the Lambda control box.
2. If you are using the serial interface, attach the 9-pin communications cable to the port labeled "serial" on the back of the control box, and the opposite end to an available com port on your computer (take note of which port you have connected it to).

Note: a 9-pin strait thru serial cable is required. If you are using the parallel interface attach the 25-pin parallel cable to the port labeled parallel (LPT) on the controller and the opposite end to an available LPT port on your computer (take note of which port you have connected it to).

If you are using USB you must install the USB driver supplied by Sutter instruments then attach your USB cable to the USB port on the Lambda controller and the opposite end to an available USB port on your computer.

3. Attach the 15-pin communication cable supplied with your Lambda controller to the connector labeled "Wheel" on the back of the control box.
4. Connect the opposite end of the 15-pin cable to the filter wheel and shutter unit.
5. Connect the power cord to an appropriate outlet and power on your Lambda unit using the power switch located on the back of the control box.
6. Switch the unit to local mode by pressing the local key located on the front of the control box (if you do not have a "local key" refer to your owners manual).
7. Cycle the wheel and shutter by pressing the "shutter" key on the key pad or use the numeric keys labeled 1 thru 9 to rotate the wheel to the correlating numeric wheel position.
8. If the unit is working properly, switch the unit to online mode by pressing the online key located on the front of the control box.

CONFIGURING FOR USE WITH YOUR META IMAGING SERIES SOFTWARE

Complete the following procedure to load the Meta Imaging Series device driver and configure the Lambda controller unit.

1. From the Windows desktop, Click Start>Programs>Meta Imaging Series> Meta Imaging Series Administrator. The Meta Imaging Series System Administrator opens.
2. Click Configure Hardware. The Configure Hardware dialog box opens.
3. Click Install System Devices. The Install Hardware dialog box opens.
4. Select Lambda 10\10-2\10-3 from the Available Hardware list and click Install. The Lambda 10\10-2\10-3 will appear in the Installed Devices list.

ARTICLE #
T20037

PRODUCTS
MetaMorph® 6.0 to 6.3

CREATED
4-Mar-2005

LAST UPDATED
4-Mar-2005

NOTES

The Lambda 10\10-2/103 driver is the correct driver for the all Sutter filter and shutter units.

5. Click Lambda 10\10-2/10-3 in the Installed Devices list to select it and click Settings. The Lambda 10/10-2/10-3 Settings dialog box opens.
6. Select your Lambda type from the Controller Type field.
7. Ensure that the appropriate filter wheels and shutters are checked (only components attached to your controller should be checked).
8. Select the port type your Lambda is connected to from the Communications Port drop-down menu.
9. Click *OK* to close the Lambda 10/10-2/10-3 Settings dialog box.
10. Click *Apply*, then click *OK* to close the Install System Devices dialog box.
11. In the Configure Hardware dialog box, click *Configure Devices*. The User Settings dialog box opens.
12. Select Lambda 10/10-2/10-3 from the Available Devices list and click *Add>>* to add it to the Claimed Devices list.
13. Select each filter wheel and label the filter positions by clicking on settings.
14. Adjust the shutter delay by selecting the shutter and clicking settings.
15. Click *Apply*, then click *OK* to close the User Settings dialog box.
16. Click *OK* to close the Configure Hardware dialog box, then Click *OK* to close the Meta Imaging Series Administrator.

© 2005 Molecular Devices Corporation