

Installing and Configuring QImaging IEEE-1394 Cameras

Abstract

Meta Imaging Series 5.0, or later, supports QImaging cameras that interface to the computer via IEEE-1394. This includes the Retiga EX, Retiga 1300, QICam, and the MicroPublisher.

Instructions

Please follow these steps to install & configure the camera:

1. DO NOT CONNECT THE CAMERA YET!
2. Ensure that your computer's VGA display settings are set for 32-bit color. This is important if you will be imaging in color. This can be done by using the Display Control Panel.
3. Shutdown the computer. Install the 1394 FireWire card supplied with your QImaging camera into an open PCI slot. Restart the computer.
4. Upon rebooting, Windows 98 Second Edition, Windows 2000, or Windows XP will automatically install the appropriate drivers. Start the Device Manager in the System Control Panel, and navigate to the "1394 Bus Controller" – the status there will indicate whether the card is working properly.
5. If your camera came with QImaging software, insert that CD now. If not, you can run the "QCapture Suite Installer" that is in the QImaging folder on the Meta Imaging Series 5.0 CD-ROM. If neither of those are available, the QCapture Suite can be downloaded from <http://www.qimaging.com/support/downloads/index.php?cat=software&subcat=windows>.
6. Once the installation screen appears, click on "Qcapture2" to begin installation
7. A setup window will open. Follow the installation instructions and choose the default choices by clicking "Next"
8. Connect the QImaging camera to the IEEE 1394 (FireWire) PCI card (either port will work) with the included IEEE 1394 cable and turn the camera on.
9. Once the installation has completed, run the program "QCapture" that is in the "QCapture Suite" menu which can be found on the "Programs" menu of the START button. The QCapture program lets you verify the camera is operational. When done, exit the QCapture program.
10. Start the Meta Imaging Series Administrator. Press Configure Hardware and then press Configure Acquisitions. Under the Available Drivers list, add the QImaging Camera Driver.
11. Click "Configure" when the QImaging Camera Driver is selected.
12. The QImaging Driver Dialog will appear.
13. Enter an identifying name for your camera or cameras in the Camera Information Tab.
14. Check "Camera has a color filter wheel" if applicable.
15. Click "OK" out of the QImaging Camera Driver Dialog.
16. Click "OK" out of the Configure Acquisition Dialog.
17. Click "OK" out of the Configure Hardware Dialog.
18. Click "OK" out of the Meta Imaging Series Administrator.

Document ID
T20010

Product
MetaMorph
MetaFluor
MetaVue
Discovery-1

Platform
Windows 95/98
Windows 2000

Created
17-May-2002

Last Reviewed
26-Mar-2003

Keywords: QImaging, Retiga EX, Retiga 1300, QICam, MicroPublisher, IEEE-1394, Firewire
Issue Type: Hardware

19. Installation of the QImaging Series Camera is complete.

Notes:

- We recommend that you use the IEEE-1394 card that came with the camera. We do not recommend using a generic or 3rd party IEEE-1394 card.
- If you install the QImaging 1394 card with the device plugged in before installing the Qcapture software, a popup icon will appear in the Taskbar prompting you to Unplug or Eject Hardware.
- After the QCapture driver is installed, a Microsoft Digital Signature Dialog will appear prompting you to continue the install or cancel. Click YES to continue, and then finish the install process.

References

QImaging may have additional installation notes at <http://www.qimaging.com/>