

QImaging camera installation

Overview

- Our QImaging driver supports QImaging cameras that use FireWire interfaces. We install [QCapture Suite](#) software to provide the camera's Windows-level driver, essential libraries, and the *QCapture* confidence test.
- A given version of *Meta Imaging Series* software must be paired with a specific version of *QCapture Suite*.

ARTICLE #
T20068

PRODUCTS
MetaMorph®
MetaFluor®
MetaVue®

CREATED
29-April-2008

LAST UPDATED
29-April-2008

<i>Meta Imaging Series</i> Version ...	requires <i>QCam Driver</i> Version ...	which is provided with ...
6.1R0 through 6.1R3	1.60	<i>QCapture Suite</i> Version 2.60
6.1R4	1.66	<i>QCapture Suite</i> Version 2.66
6.1R5 through 6.2R0	1.68.2	<i>QCapture Suite</i> Version 2.68.2
6.2R1 through 6.2R2	1.68.4	<i>QCapture Suite</i> Version 2.68.4
6.2R3 through 6.3R0	1.68.6	<i>QCapture Suite</i> Version 2.68.6
6.3R1 through 6.3R2	1.71	<i>QCapture Suite</i> Version 2.71
6.3R3 through 6.3R7	1.73	<i>QCapture Suite</i> Version 2.73 and <i>QCapture Pro</i> 5.1.1.14 .
7.0R0 through 7.0R3	1.81	<i>QCapture Suite</i> 2.81
7.0R4 through 7.1.7	1.90	<i>QCapture Suite</i> 2.90
7.5.0 through 7.5.3 32-bit	2.05	<i>QCapture Suite</i> 2.95
7.5.0 through 7.5.3 64-bit	x64-2.05	<i>QCapture Suite</i> x64 2.95

- Do not attach more than one QImaging camera to a given FireWire adapter.

Installation instructions

1. Install the FireWire card from QImaging;
2. Install *QCapture Suite*;
3. If the camera uses an external power supply, turn the camera on;
4. Attach both ends of the FireWire cable that runs between the camera and the FireWire port; this allows the FireWire hardware to handle its part of the installation;
5. The *Found New Hardware* Wizard starts to allow Windows to configure itself to see the camera. Allow the Wizard install software automatically.
 - a. (*Windows XP Professional with Service Pack 2*) If the Wizard offers to check Windows Update for the most recent driver, choose "Not this time".
6. Configure *Meta Imaging Series* to acquire from the camera.

Upgrading QCapture Suite for an existing camera

1. Make sure the camera is powered up;
2. Right-click the *Windows Device Manager* entry for the camera, which brings up a pop-up menu containing *Update Driver*; *Disable*; *Uninstall*; *Scan for hardware changes*; *Properties*.
3. Choose *Uninstall* and confirm that you really wish to do that. This notifies Windows that the camera is to be removed and information about the camera excised;
4. Unplug either end of the cable that connects the camera to the FireWire port. This allows the FireWire hardware to handle its part of the removal of the hardware.
5. Use the *Add or Remove Programs* control panel to uninstall the QImaging software:
 - a. Uninstall the imaging application. It might be listed as *QCapture Suite* or *QCapture Pro*.
 - b. Uninstall any QImaging camera driver.
6. Install the required version of *QCapture Suite*.
7. If the camera has its own power supply, make sure the camera is powered up;
8. Plug the end of the cable back in; this allows the FireWire hardware to handle its part of the installation, and triggers the *Found New Hardware* Wizard;
 - a. (*Windows XP Professional with Service Pack 2*) If the Wizard offers to check Windows Update for the most recent driver, choose "Not this time".
9. Configure *Meta Imaging Series* to acquire from the camera.

If there are any further questions, please contact Technical support at 800-635-5577 x1820 or support.dtn@moldev.com.