

Photometrics Camera Troubleshooting

Document ID
T10245

Created
17-Aug-2000

Abstract

Following are some recent problems, and some ways we have found to work around them.

Instructions

Problem: Meta Imaging Series software "locks up" when started, if the camera is off.

On some systems, Meta Imaging Series applications will "lock up" at startup if they are configured to use a Photometrics camera, and the Photometrics camera is not on when the Meta Imaging Series application tries to start up.

Solution

The "lock up" problem occurs if the PVCAM.INI file specifies more than one camera. One way to solve this problem is to use Find Files and search for all copies of PVCAM.INI on the startup drive (usually C:) and then delete all of those files. After doing so, run the PCSETUP95 (win95/98 systems) or PCSETUPNT (winNT systems) that is in the Photometrics program group. Re-enter the name of your camera. Run Video Manager and reconfigure the Photometrics driver.

Once those steps have been taken, Meta Imaging Series applications will be able to determine if the Photometrics camera is not powered on at startup, and they will be able to put up a message asking you to power on the camera before continuing.

Problem: Camera error 53 at startup.

Camera error "53" at startup. Camera will not work with Photometrics CAMTEST application.

Solution

If you are using a PCI Model 2 board: The original PXL camera and the new Quantix camera have the same type of connector, but use a different cable to interface to the PCI Model 2 board. Not only do they use a different cable, but the cable is direction-specific. Please verify with Photometrics that you have the correct cable. If it does not work in one orientation, then swap the cable and retry.

Keywords: photometrics hardware
Issue Type: Hardware