

Mutech MV-1500 Configuration Notes

ABSTRACT

The MV-1500 is a 16bit digital-only frame grabber from Mutech. It is intended for customers who only have digital cameras. If control of both digital and video cameras is desired, the recommended solution is the installation of both a MV-1500 and a FlashBus card. The MV1500 is typically used to communicate with Hamamatsu digital cameras.

CABLE NOTES

- There is one digital cable that supports all Hamamatsu cameras.
- The cable can be provided by Molecular Devices Corporation or ordered from Cables To Go as part # CCUNIVMV-1500.
- The cable is **unidirectional** so be sure to install it in the proper direction. One end is labeled "Camera" and the other "MV1500". If these are connected in reverse then the camera will not work with Meta Series software.

DRIVERS

When the MV-1500 board is first installed, Windows will detect it. When Windows prompts for a driver, point it to the "MuTech MV1500\win2k" folder on the Meta Imaging Series CD-ROM. This folder contains the proper driver files for the board installation.

ADMINISTRATOR ACCESS

The MV-1500 driver requires Windows login Administrator rights to function correctly. If Administrator rights are not available then the driver will be unable to allocate memory necessary to stream frames or view fast live images.

If you must login to Windows using an account other than Administrator, then the following procedure must be used to enable proper access for the other accounts:

- Authorized user should login to Windows using the Administrator account.
- Open the control panel called Administrative Tools
- Launch the program within Administrative Tools called Local Security Policy
- Open the Local Policies folder
- Open the User Rights Assignment folder
- Highlight the Increase Scheduling Priority entry and double-click it
- Click the Add Users or Group button
- Enter each user account which needs to run the MV-1500 driver
- Click OK to complete the procedure

ARTICLE #
T10168

PRODUCTS
MetaMorph® 4.6 or later
MetaFluor® 4.6 or later

CREATED
18-Dec-1998

LAST UPDATED
04-Sep-2005