

## Flashbus Driver for Windows 2000

**Document ID**  
T10250

**Created**  
03-Jul-2001

### Abstract

The Flashbus driver has been updated to work with Windows 2000. To use the updated driver, you must have Meta Imaging Series version 4.6 with the 4.6.4 (or later) update installed. The 4.6.4 update is the first update to contain the fully Windows 2000 compatible driver. There are two different installation routines. One applies if you are using the Flashbus for the first time under Windows 2000, while the other applies if the board is already installed in the Windows 2000 computer.

### Instructions

#### Using the Flashbus for the First Time under Windows 2000

- ◆ Do not install the board
- ◆ Install Meta Imaging Series 4.6
- ◆ Download and install the latest 4.6 update, currently 4.6.4. You can find the update by clicking on this link: [T10249](#).
- ◆ Shutdown the computer.
- ◆ Verify that the Flashbus board is jumpered correctly. You can find the jumper information by clicking on this link: T10165.
- ◆ Install the Flashbus board in the computer.
- ◆ Startup the computer. Windows 2000 will detect the Flashbus board. Identify the board to Windows by selecting the "fbusmv2k.inf" file that is located in the \mm\input\flashbus\Win2K\ directory.
- ◆ Once Windows 2000 has recognized the board, you can configure the board using the Administrator.
- ◆ Note that after you configure the board, you will be prompted to reboot the computer. This allows the driver file to get installed into the c:\winnt\system32\drivers directory.

#### Using the Flashbus if the Board is Already Installed in a Windows 2000 PC

- ◆ Install Meta Imaging Series 4.6
- ◆ Download and install the latest 4.6 update, currently 4.6.4. You can find the update by clicking on this link: [T10249](#).
- ◆ Open the Control Panel and start the System control panel.
- ◆ Choose the Hardware tab, then press the Device Manager button.
- ◆ Under "Sound, video and game controllers" you will find the "FlashBus MV" driver entry.
- ◆ Double-click on the "FlashBus MV" driver and choose the Driver tab. Press the Update Driver button.
- ◆ Using the Upgrade Device Driver Wizard, search for the "fbusmv2k.inf" file that is located in the \mm\input\flashbus\Win2K\ directory.
- ◆ After Windows has updated the driver, run the Meta Imaging Series Administrator.

**Keywords:** flashbus hardware  
**Issue Type:** Hardware

- ◆ Reconfigure the Flashbus board, and then reboot the computer.