

Meta Imaging Series Online Support

[Home](#) · [Software Updates](#) · [Supported Hardware](#) · [Technical Notes](#) · [Contact Us](#)

→ [Technical Notes](#) → [Browse Technical Notes](#)

Use the **small arrows** to the left and right of the column name to **sort** this list in ascending or descending order.

↑ ID ↓	↑ Title ↓
T20050	Meta Imaging Series information and support resources
T20049	Creating an 8-bit copy of 16-bit images in MetaMorph
T20048	An introduction to MetaFluor
T20047	MetaFluor digital triggers
T20046	Using the TILL Photonics Polychrome V with Meta Series 6.3r3 and beyond
T20045	Using the Multi Wavelength Cell Scoring application module in MetaMorph
T20044	Creating custom toolbars, menu and icons in MetaMorph and MetaXpress
T20043	Using the Angiogenesis Tube Formation application module in MetaMorph version 6.3 and higher
T20042	Using the Count Nuclei application module in MetaMorph version 6.3 and higher
T20041	Zeiss AxioImager.Z1 microscope support in version 6.3r2 and later
T20040	Using the Morphology Filters
T20039	Spatially calibrating images in MetaMorph
T20038	Installing Diagnostic Instruments SPOT cameras with MetaMorph Versions 6.0 through 6.3
T20037	Installing and configuring Sutter Instruments Lambda filter wheel/shutter devices
T20036	MetaMorph application modules overview
T20035	How to use image file variables to save images in a journal without user interaction
T20034	How to Use Variables to Create Reusable Screening Analysis Journals for any Wavelength Combination
T20033	Advanced Journaling Tools Available on the Actions Tab of the MetaMorph Journal Editor
T20032	Using the Universal Imaging Corporation FTP Site
T20031	Correcting for Uneven Background Using Morphology Filters
T20030	Using Variables to Override Function Settings in a Journal
T20029	Running the Neurite Outgrowth Dialog Box in MetaMorph Premier
T20028	Eliminating Memory Key Driver Conflicts with Meta Imaging Series Hardware
T20027	PIFOC E-662 RS-232 Serial Cable
T20026	Important Differences Between Previous Versions of 6.1 and Version 6.1r6 and Above
T20025	Using Integrated Morphometry Analysis (IMA) to select parts of an image for further analysis
T20024	Custom Data Logging in MetaMorph
T20023	Using the KeyPad in Meta Imaging Series Applications
T20021	Using AutoQuant Deconvolution
T20020	Microsoft SQL Server 2000 Desktop Engine (MSDE 2000): Installation Notes, Configuration Notes, and Issues for the Image B...
T20019	Zeiss Automated Microscopes and the Light Manager
T20018	Installing Microsoft SQL Server Desktop Engine Service Pack 3
T20017	Improving Signal To Noise When Imaging Fura2 with a Monochromator
T20016	Using the Measured PSF 3D Deconvolution Command
T20015	Troubleshooting the Image Browser
T20014	AttoFluor AIC 4-Position Filter Changer
T20013	MetaMorph Functions whose Parameters can be Overridden by Variables
T20012	Polychrome IV Configuration for MetaMorph 5.0
T20011	Hamamatsu C4880-81 Horizontal Offset
T20010	Installing and Configuring QImaging IEEE-1394 Cameras
T20008	EFOS/EXFO X-Cite
T20006	Multi Dimensional File Format
T20005	Prairie Technologies AOTF Driver
T20004	Nikon Eclipse TE-2000 Configuration
T20003	Hamamatsu Orca Cable Information
T20002	Screen Acquisition option is not present
T20001	The use of journals and variables during Multi Dimensional Acquisition
T10258	Diagnostic Instruments SPOT RT 3.4 Driver Update
T10257	Photometrics PVCAM 2.5.8 Driver Update
T10256	Using the CAIRN Optoscan with DAC Interface
T10255	Using the ComputerBoards PCI-DDA02 with Windows 2000
T10254	Linkam Temperature Controlled XY Stage
T10253	Using the IO Tech DaqBoard/2000 Device
T10252	Installing and Configuring Hamamatsu IEEE-1394 Cameras
T10251	MV-1500 Driver for Windows 2000

[T10250](#) Flashbus Driver for Windows 2000

[T10249](#) Cairn Optoscan Cabling Information

[T10248](#) Ludl Serial Cable

[T10245](#) Photometrics Camera Troubleshooting

[T10244](#) Photometrics NORMAL vs. ALT Modes

[T10243](#) MetaMorph STK and TIFF File Formats

[T10240](#) SPOT Camera - Uninstalling and Reinstalling the SPOT Software

[T10239](#) Windows [VCACHE] - Saving Images "eats up" RAM

[T10238](#) Prior XYZ - Images Acquired Before Motor Stops

[T10237](#) Quantix 57 Usage Notes

[T10236](#) Installing the CRI VariSpec RGB or the CRI MicroColor

[T10232](#) Tools for Determining Computer Resource Allocation

[T10231](#) Converting Spot Camera images to the correct greyscale range

[T10230](#) Adobe PhotoShop can't open 16-bit TIFF images, or displays them as black

[T10229](#) 8x Binning with the Hamamatsu ORCA Camera

[T10228](#) SPOT RT Camera

[T10226](#) Memory Key and Parallel Ports -- Preventing Conflicts

[T10224](#) Problems with auto-exposure and Photometrics cameras

[T10223](#) Meta Imaging Series Updater: 4.0 to 4.1.7

[T10222](#) Nikon Remote Focus Accessory - Configuration Information

[T10220](#) Data Logging to MicroCal Origin

[T10218](#) Photoshop v.5 User Caution

[T10211](#) Dialog Box Appearance

[T10208](#) Olympus AX-70 Driver Update and Configuration Information

[T10200](#) Using Classifiers to Exclude Unwanted Objects From Colocalization Measurements

[T10199](#) Measuring or counting objects within objects

[T10198](#) Counting and Measuring Branch Points on Tree Like Structures

[T10197](#) Using Variables in MetaMorph to Create Concentric Regions

[T10196](#) Sutter Lambda 10, Lambda 10-2 Device Not Responding Error Message

[T10191](#) Installation Issues for Meta Imaging Series 4.0

[T10190](#) FlashBus Trigger Out and Switch Input

[T10188](#) Display Option: Overlay Image Command

[T10187](#) PI Driver Update for Meta Imaging Series 4.0

[T10185](#) Meta Imaging Series - Windows Compatibility

[T10184](#) Ludl Driver - Advanced Features

[T10182](#) Lambda Utility to Initialize / Test the Lambda 10-2

[T10181](#) Zeiss MTB Microscope

[T10180](#) Using the Princeton Instruments CCD57 Back Illuminated Frame Transfer CCD Camera with Meta Imaging Series 3.6

[T10179](#) Meta Imaging Series Region File Format

[T10178](#) Using the Roper Scientific COOLSNAP with MetaMorph Version 4.0 and Above

[T10177](#) The Sutter Lambda-10DE Fast Filter Wheel

[T10176](#) Using the Orca II with Meta Imaging Series 4.0 and Above

[T10170](#) Nikon E1000 Serial Communications

[T10168](#) Mutech MV-1500 Configuration Notes

[T10167](#) Updated MuTech Driver Files for Meta Imaging Series Version 3.6

[T10166](#) PI VInput Driver for Meta Imaging Series Version 3.6

[T10165](#) Using the FlashBus Card with Meta Imaging Series 3.6 and 4.0

[T10164](#) MuTech Analog Settings

[T10163](#) Uniblitz Shutter Serial Cable

[T10161](#) PI Interline Temperature Settings

[T10160](#) Using Photometrics Cameras with Meta Image Series Versions 3.5 and Above

[T10158](#) Orca Exposure Times: Streaming vs. Single Frame Acquisition

[T10157](#) Configuring a MetaMorph System to Run as an Acquisition or an Offline System

[T10156](#) Adaptec Easy CD Creator Software Problems Using Teac CDR-55S Drives

[T10155](#) Adjusting the Horizontal Offset for Hamamatsu C4880 Cameras

[T10151](#) Installing Meta Imaging Series 3.6 on Windows NT Systems

[T10150](#) 1-2-3 97 Does Not Accept Logging by Meta-Series Products

[T10146](#) DDE to Microsoft Excel 97 crashes Excel

[T10144](#) Updated FlashBus Driver Files for Meta Imaging Series 3.6

[T10143](#) User Call command (for Visual Basic programs) won't run

[T10140](#) MuTech Device Manager File

[T10135](#) Updated Diagnostic Instruments SPOT Driver for MetaMorph 3.5

[T10133](#) Measuring the extension and retraction of cells or the movement of cells using Measure Colocalization

[T10132](#) PI VInput Driver for MetaMorph 3.5

[T10131](#) Anti-virus Program Shows ".OUT" Files Infected

[T10128](#) The Format of IMG Files

[T10127](#) Using the PolyChrome Monochromator to Shutter Light

[T10126](#) How to Determine the Version of Windows 95 that Your Computer Uses

[T10125](#) Using classifiers to select objects for measuring brightness over time.

[T10124](#) Meta Series Software Parallel Port Pin Numbering

[T10122](#) Ludl MAC 2000 / MAC 5000 Controller Setup

[T10121](#) Updated MuTech MV-1000 Driver Files for MetaMorph 3.5

[T10120](#) ORCA-100 produces black images when focusing

[T10119](#) Imena PC90 SCSI Adapter Switch Settings

[T10118](#) Iomega PC2 SCSI Adapter Switch and Jumper Settings

[T10115](#) Integrating Camera Q&A

[T10114](#) Updated Photometrics Driver Files for Meta Imaging Series 3.5 and 3.6

[T10113](#) NI-DAQ Software Updates from National Instruments

[T10108](#) Video Output from the ATI 3DPRO TURBO PC2TV Card

[T10107](#) Using The Multi-User Chooser with Version 3.5 of Meta-Series Software

[T10106](#) Leica DM RXE and IRBE Cable

[T10104](#) Universal Imaging SCSI FAQ

[T10099](#) PTI DeltaRAM Installation & Testing Notes

[T10098](#) Hardware Integration Enhancement Option for MuTech

[T10097](#) Physik Instruments PIFOC Z Motor Interface Hardware

[T10095](#) Polychrome Monochromator Interface Hardware

[T10094](#) MuTech MV-1000 Troubleshooting Q&A

[T10093](#) Installing the Automated Microscope Option in

[T10092](#) Errors Using Protected Zip Disks on a Computer with an Adaptec SCSI Controller

[T10089](#) Workaround for Zip Tools Conflict with MetaMorph, MetaFluor

[T10088](#) Equipment Needed to Install a Zip Drive on a Universal Imaging System

[T10086](#) The Iomega Zip Drive and UIC Products: Introduction and Profile

[T10082](#) MuTech Jumper Configuration for Hamamatsu Digital Camera

[T10079](#) Transfer RGB image to video using Matrox MVP board

[T10078](#) Common Problems With Ratio Imaging Experiments

[T10077](#) Using Classifiers to Exclude (or Include) Unwanted Objects From Region Measurements

[T10076](#) Hamamatsu C4880-80 Delay When Binning or ROI Changed

[T10072](#) How to install a Photometrics Camera for use with Meta Series Software 3.0

[T10070](#) Configuring the MuTech MV-1000 Video Acquisition Board Driver in Video Manager

[T10069](#) Using the MuTech MV-1000 with the Hamamatsu 4742-95 or C4880-80/81

[T10068](#) Using Digital and RGB cameras with the MuTech Board

[T10067](#) Using the MuTech Board with MetaFluor or MetaGFP

[T10066](#) MuTech Camera Configuration Files

[T10065](#) MuTech Cable Specifications

[T10064](#) MuTech Hardware Configuration Comparison

[T10063](#) MuTech Board Jumper Configuration for Analog Acquisition

[T10062](#) Matrox Millenium 16-bit Driver for Windows 95

[T10058](#) Adaptec SCSI Controller Driver Updates

[T10057](#) Using the Pinnacle RCD-5040 or RCD 4x4 with an Adaptec 2910 or 2940 Controller

[T10056](#) Using MetaMorph for FISH and CGH

[T10055](#) Marzhauser Stage Connection Test Procedure

[T10054](#) How is "Standard Area" used to determine a count?

[T10053](#) Installing Hewlett-Packard LaserJet disables MetaMorph

[T10051](#) Marzhauser MultiControl 2000 Cable Information

[T10050](#) Ludl Stage Motors Not Reporting Their Positions

[T10049](#) Logging Data to Non-English Versions of Excel

[T10048](#) Diagnosing Uniblitz Shutter Problems

[T10047](#) Wait for Digital I/O Accuracy

[T10046](#) Sutter DG4 Control Error

[T10043](#) Hamamatsu C5985 Firmware

[T10042](#) Windows 95 and Windows NT 4.0 support for Polaroid CI3000S, CI5000 Digital Palette devices

[T10041](#) Configuring Image-1 or MetaMorph to control the LQ4000 OMDR

[T10038](#) Computer System Driver and BIOS Updates

[T10036](#) AT200 Control of an External Shutter

[T10033](#) Orca and Orca 2 Gain Settings

[T10032](#) INI command failed

[T10031](#) Installing 3COM 3C905/3C905B Fast EtherLink XL Adapters

[T10030](#) Uniblitz Shutter Cables

[T10029](#) Windows Hardware Compatibility List

[T10028](#) Matrox MGA Video Driver Updates

[T10027](#) Adaptec I/O Interactive Database

[T10022](#) Using MetaMorph to align wavelengths in MetaFluor or MetaGFP

[T10021](#) Viewing "Live Video" with MetaFluor Version 3.5 and Above

[T10020](#) Configuring Windows 95 for the Image-LC

[T10019](#) Configuring Windows 95 for the MVP/AT

[T10017](#) An Introduction to MetaFluor

[T10016](#) MetaFluor Quick Titration Calibration

[T10015](#) Using MetaFluor to Analyze Experiments Stored on OMDR

[T10014](#) MetaFluor and the MVP Video Driver

[T10013](#) Aligning Mosaic Montages

[T10012](#) Using Optical Density (Scaled) for Densitometry

[T10011](#) Meta Configuration Files

[T10009](#) Shading Correction and Background Subtraction

[T10005](#) Analog measurements option for MetaFluor

[T10004](#) Maximum number of gray levels displayed under Windows

[T10003](#) MetaFluor IMD display mode

[T10002](#) MetaFluor Digital Triggers

[T09014](#) Updated Matrox MVP-AT Driver Files for MetaMorph 3.5

[T09012](#) Photometrics Driver Files for MetaMorph 3.0
[T09011](#) PI VInput Driver for MetaMorph 3.0
[T09010](#) Matrox MVP-AT VInput Driver for MetaMorph 3.0
[T09009](#) MuTech MV-1000 Driver for MetaMorph 3.0
[T09002](#) Meteor VInput Driver for MetaMorph 3.0

[Top](#)