

Transferring Meta Imaging Series Installation across multiple PCs

Overview

- Pre-installation requirements
- Original system Meta Imaging Series configuration data
- Clean installation of Meta Imaging Series on the target computer
- Transfer configuration data to target computer
- Make final edits to configuration data on target computer

ARTICLE #
T20097

PRODUCTS
MetaMorph®
MetaFluor®
MetaVue®

CREATED
11-Aug-2008

LAST UPDATED
20-Aug-2008

Pre-Installation Requirements

This document is for transferring configuration data from one Meta Imaging Series installation to another. The requirements for this transfer are listed below:

- The transfer is from the same version of the Meta Imaging Series software (e.g. 7.5 to 7.5)
- The transfer is from the same System ID of the Meta Imaging Series software (e.g. AAAAA to AAAAA). This document will not be accurate for attempting to transfer configuration data from one System ID to a different System ID (e.g. AAAAA to BBBB).
- All third party software and hardware must be installed and recognized by the target system - where applicable - before the transfer is attempted. (e.g. third-party SDK, control software, PCI cards and drivers, etc.) For additional questions, see hardware specific technical notes.

Original system Meta Imaging Series configuration data

Key configuration data for Meta Imaging Series software includes the following:

Meta Imaging Software configuration data consists of the following files and folders:

UICSHELL.INI – contains license information about the installation

MMADMIN.INI – contains user / password information

MM\Groups\ folder – contains individual user configuration data

MM\Hardware\ folder – contains system hardware configuration data

MetaMorph.INI, MetaFluor.INI, MetaVue.INI, etc. – application group specific information. The installation will have at least one of these files, possibly more depending on which applications are available.

In addition, there are user-defined configurations that must be moved as well, such as:

Illumination settings (.ILS)

Magnification settings (.MGS)

Acquisition State settings (.ACQ)

Calibrations (.CAL)

Saved Regions (.RGN)

ND State files (.ND)

Taskbars (.JTB)

Journals (.JNL)

All of the above user-defined settings, except for the journals, are stored in temporary configuration (.cfg) files within the individual user's folder in the main MM\Groups folder by default. If those configuration files are deleted then the user settings are lost. If these files are present, then copying the MM\Groups\ folder is sufficient to move the user settings from one installation to another¹.

Any Journals created must be copied from the original system to the corresponding locations within the clean installation. Failure to do so may cause the journals to malfunction.

Clean installation of Meta Imaging Series software

Install a clean version of the Meta Imaging Series software into a home directory that is identical to the original system. That will ensure the least amount of disruption and re-configuration needed.

Transfer configuration data to target computer

Once the new PC hardware has been installed and verified, copy the above files and folders to the corresponding locations in the clean installation, overwriting the existing folders and files.

Make final edits to configuration data on target computer

Open the Meta Series Administrator and click on Configure Hardware. Review all hardware configured and make any changes necessary, including communication ports.

After verifying the hardware in the Meta Series Administrator, open MetaMorph and review the user settings listed above. Make any changes necessary.

¹ To prevent this loss of data, the user can store their settings to the hard drive as backup files. (These files are labeled with the file extensions listed above.) To create a backup file, the user must click on the Backup button from whatever settings window they are saving. For example, if the user wants to create a backup of the Illumination Settings, then they would click on the Backup button in the Configure Illumination Settings window.

Once all of the settings have been backed up, the user can then copy the files from the existing installation and paste them in the corresponding location in the new installation. To restore the settings, select Restore from the settings window in the new installation.