

Converting a Meta Imaging Series® software Installation from one Installed system ID to another system ID of similar options

Overview

- Pre-Transfer requirements
- Clean installation of Meta Imaging Series® software on the target computer
- System should be configured and up and running
- Backing up files and converting system
- Finalizing your conversion

Pre-Transfer Requirements

This document is for modifying a Meta Imaging Series® software installation so that it can be run with another dongle containing a different system ID

- Functional installation of your Meta Imaging Series® software configured and able to control all hardware available.
- The transfer is only possible for system ID's that are able to run the same versions of the Meta Imaging Series® software (e.g. 7.6 to 7.6)
- The transfer should only be attempted with system ID's that are of the same category of Meta Imaging Series® software (e.g. Premier to Premier) or (e.g. Basic to Basic)
- Registration code for new system ID and Authorization Code for new system ID

Clean Installation of Meta Imaging Series® software

Install a clean version of the Meta Imaging Series® software into any location of your choice. Please note the root path of your installation for future reference.

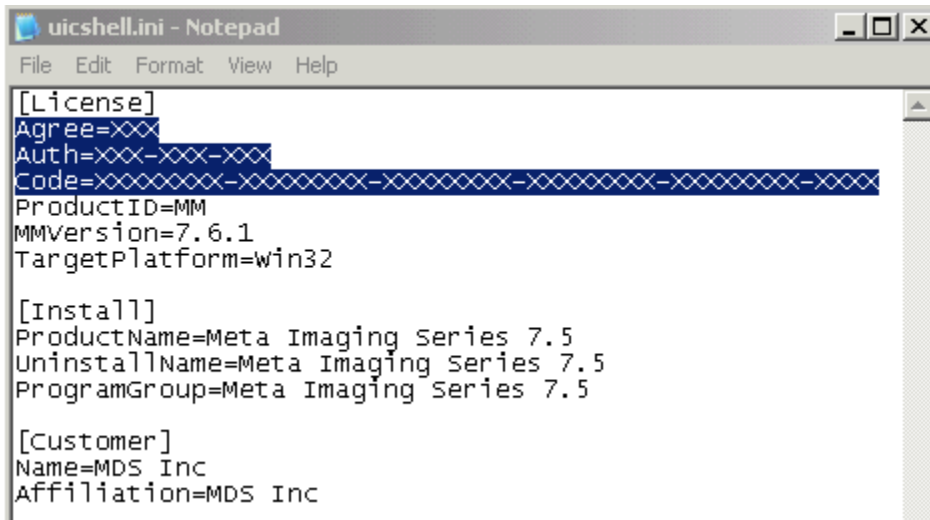
System should be configured and up and running

Complete any available updates to the Meta Imaging Series® software installation using the Meta Imaging Series® software Updater. Install all required 3rd party drivers for your hardware and plug-in and configure your hardware. Make sure that all hardware is functional.

Backing up files and converting system ID's

- Go to the root directory of your Meta Imaging Series® software installation and find the UICshell.ini file.
- Make a backup copy of this file into a safe location.
- Open the original UICshell.ini found in the Meta Imaging Series® software root directory using Notepad.exe for editing.
- Edit the following lines in the file with the updated information fro your new system ID
 - (e.g. Agree=XXX to Agree-123)

- (e.g. Auth=XXX-XXX-XXX to Auth=123-456-789)
- (e.g. Code= XXXXXXXX-XXXXXXXX-XXXXXXXX-XXXXXXXX-XXXXXXXX-XXXX to Code=12345678-12345678-12345678-12345678-1234)



```
uicshell.ini - Notepad
File Edit Format View Help

[License]
Agree=xxx
Auth=xxx-xxx-xxx
Code=xxxxxxxx-xxxxxxxx-xxxxxxxx-xxxxxxxx-xxxxxxxx-xxxx
ProductID=MM
MMVersion=7.6.1
TargetPlatform=win32

[Install]
ProductName=Meta Imaging Series 7.5
UninstallName=Meta Imaging Series 7.5
ProgramGroup=Meta Imaging Series 7.5

[Customer]
Name=MDS Inc
Affiliation=MDS Inc
```

Finalizing your conversion

- Unplug your old dongle and insert the new one containing the new system ID
- Now run the Meta Imaging Series[®] software Updater using the last update that was completed as the source file during your update. Please make sure that the correct system ID is listed during your update.
- Your system should now be setup for the new dongle. Please retest all hardware and options that are available for the new system ID. Only options for the new system ID will now be available.