

Zeiss Axiomager.Z1 microscope support in Version 6.3r2 and later

ABSTRACT

Meta Imaging Series 6.3r2 and up now supports the Zeiss Axiomager.Z1 automated microscope. Support for this microscope requires installation and configuration of the Zeiss **MTB 2004** microscope software toolkit.

REQUIREMENTS

- Meta Imaging Series 6.3r2 or later.
- Appropriate device control options.
- Zeiss Axiomager.Z1 microscope.

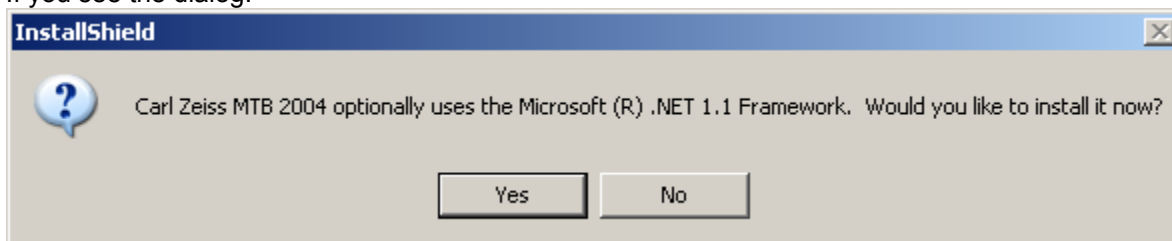
OVERVIEW

The first step to controlling the Zeiss Axiomager.Z1 microscope from MetaMorph is to install the Zeiss MTB 2004 toolkit. This toolkit can be found on our website at <http://www.universal-imaging.com/ftp/support/hardware/ZeissMTB2004/Ver1021/>.

There are two zip files available from the FTP link. If your computer already has Microsoft .NET frameworks 1.1 installed, then you only need to go to download the "MTB2004-RDK1_0_2_1wo_net_Runtime.zip" file, which is less than 4 MB. However, if your computer does not have .NET then you will have to download the much larger (about 26 MB) file called "MTB2004-RDK1_0_2_1.zip". We suggest you download the smaller one, try to install it, and if it asks for .NET then cancel the install and re-download the larger file and install that instead.

Once you have downloaded the zip file, unzip it to a temporary directory such as c:\temp\ and then run the "setup.exe" that is contained within. This will install the Zeiss toolbox.

If you see the dialog:



This means that the .NET framework must be installed. If you only downloaded the "MTB2004-RDK1_0_2_1wo_net_Runtime.zip" file, then you must select "No" and re-download the larger file, "MTB2004-RDK1_0_2_1.zip". If, however, you are running the setup.exe from the larger file, then answer "Yes" to have the .NET framework installed.

Leave all the options as selected (except for entering your name and affiliation) and proceed with the installation.

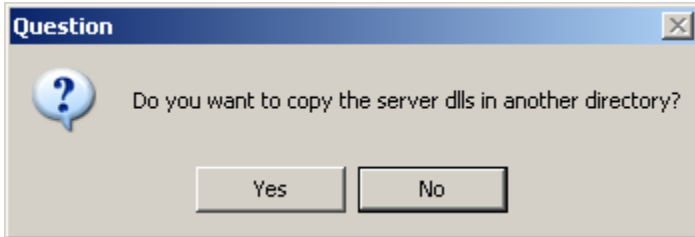
At the end of the toolbox installation, the following dialog will appear:

ARTICLE #
T20041

PRODUCTS
MetaMorph® 6.3r2 or later

CREATED
25-Jul-05

LAST UPDATED
25-Jul-05



You must answer “Yes” to this. A directory chooser window will appear hidden behind the setup screen. Find the directory chooser dialog, and choose the c:\mm directory (or wherever Meta Series is installed).

Then the Zeiss installer will ask again the same question, this time choose “No”.

Once the installer exits, use explorer and confirm that it has created the directory c:\MM\MTB 2004\. Navigate to this directory and copy the two files within (“MTBApi.dll” and “MTBApi.dll.config”) from there to the c:\MM directory.

Double-click the “MTB Configuration” icon that is available on the desktop. Use this program to configure the microscope.

Should you wish to “emulate” the microscope, then choose “New Configuration” from the file menu. Double click on it, highlight “Other Stand” and then double-click on Axio Imager from the stand choices on the right. Then click the “Simulate Hardware” checkbox that appears.

Continue through this program, configuring the proper options. Then exit the program.

Start up the Meta Imaging Series Administrator, press “Configure Hardware”, and then “Install System Devices”. Find the driver “Zeiss COM MTB” on the left and then add it to the right. This driver will detect and use all the components present in the microscope that are supported by MetaMorph.