

Updating MetaXpress[®] Software



The purpose of this document is to guide the user on updating MetaXpress software through the Meta Imaging Series Updater.

- 1. Go to the MetaXpress Updates Website
 - a. http://www.meta.moleculardevices.com/mxupdates/
- 2. Enter you System ID and click Log In
 - a. Your System ID can be found on the MetaXpress USB dongle OR
 - b. MetaXpress software in the main menu, click on **Help** > **About MetaXpress**



MetaXpress Software Updates

Welcome to the MetaXpress® Software Updates website.

This site provides update files for MetaXpress software ("Update Files"), as well as certain detailed information about those updates such as release notes. The Update Files may include error corrections, bug fixes, minor improvements, or incremental features or functionalities, but do not include any upgrades that provide substantial new features or additional functionalities ("Upgrades"). The information on this site is maintained on a regular basis to reflect the currently shipping version of the MetaXpress software.

If you are interested in obtaining a license to any Upgrade of a MetaXpress software product, please contact your <u>sales representative</u> or email us at <u>meta.admin@moldev.com</u>.

Note: Cookies and Java Script must be enabled.



To log in, enter your system ID here:

and press Log In

Log In Without ID Note: You will be unable to download updates

3. The update page with available updates for MetaXpress and Custom Module Editor will appear



- 4. Click on the appropriate link in the left hand column for the correct version of MetaXpress
- 5. You will then be redirected to a page with links for 32-bit and 64-bit versions of the software. Click on the apropriate version for your computer.
 - a. To check what version of Windows you are running, click on **Start**, right-click **Computer**, and then click **Properties**
 - b. If "x64 Edition" is listed under System Type, you are running the 64-bit version of Windows



- 6. When the license agreement appears, click **I ACCEPT THE TERMS IN THIS AGREEMENT**
- 7. The update file will be automatically downloaded to your computer



CAUTION! Do not attempt to unzip the update file. The Meta Imaging Series® Updater unzips the file during the update process

8. From the Windows Start menu, select Programs > MetaXpress 6 > Meta Imaging Series Updater



9. Click Next

10. Click on the ellipsis (...) button and the Windows Open dialog will appear

11. Browse for the update file you just downloaded

Meta Imaging Series Updater		Meta Imaging Series Updater
Select Patch File Select the patch file for this update		Select Patch File Select the patch file for this update
Patch File caelect a patch file>	_	Patch Fie C-\Users\kuciane ganiko\Dxouments\SOFTWARES_U_\ms5_3_0_5_x64_2014-11-21.zp
cBack [Next) [Cancel	(Back Next) Canal

12. Click Next

13. The Hardware Key Confirmation page will appear

Confirm that a v	ntirmation alid hardware key is attached to the system
System ID	Read System ID from Key
Key Location • Use Local Key Network Key Ser Server Address:	ver localhost

- 14. If you are using a **network key**, skip to Step 15. If you are using a MetaXpress Software hardware key to update the software on an instrument or analysis (offline) workstation computer, make sure the hardware key is inserted into a USB port, select **Use Local Key** and then click **Read System ID from Key**
- 15. If you are using a **network key** to update the software on an analysis (offline) workstation computer, select **Network Key Server**, type the server address (static IP address or full computer name) where the network key is installed, and then click **Read System ID from Key**

16. The Updater should now display your System ID in the System ID field

Note! If the software does not detect the hardware key, you will not be able to click the Next button to go the next page of the Updater. If this occurs, make sure the USB key is connected, the network key server software is installed, and that the appropriate firewall settings are applied.

- 17. Click Next
- 18. The User Information and Codes page will appear

Enter the req	uired user information and cod	es to proceed with this ins	tal		
Registered Owner	Information				
Name:	Molecular Devices				
Affiliation:	Molecular Devices				
Codes					
Registration Code:					
Authorization Code	e:				
Retrieve Codes:	Using the Internet	From a Disk	Manually		
Note:	Registration Codes and Author	orization Codes are case s	ensitive.		

19. Enter your Name and Affiliation (optional)

20. Enter the codes using one of the following methods:

- a. If the computer is connected to the internet, click **Using the Internet**. The Updater retrieves the codes and displays them in the appropriate fields.
- b. If the codes are on a disk, insert the disk and click **From a Disk**. The Updater copies the codes from the disk to the appropriate fields.
- c. If the computer is not connected to the internet, click **Manually** and type the registration and authorization codes in the appropriate fields .You can find the codes on the letter that accompanied the order or you can go to
 - i. http://www.meta.moleculardevices.com/authorize/
 - ii. You will be asked to enter your System ID.

21. Click Next

- 22. The Ready to Proceed dialog will appear. Click Update
- 23. When the installation is complete, the Update Complete screen appears

24. Click Finish

1311 Orleans Drive Sunnyvale, CA 94089 25. The Meta Imaging Series Updater opens the Meta Imaging Series Administrator screen

Group Name	Hardware Settin	g File Associatio	Select a Group and	l Pre
MetaFluor	Default			126.
MetaFluor Offline	Offline		Assign Hardwa	ſê
MetaMorph MetaMorph Ofline	Offline		Drop-ins/Toolb	ats
MetaVue	Default		Clear Setting	s
MetaXpress	Default			
•			Edit Defaults	
Pressing Set File Associati pe launched when .stk an (marked by *)	on will set the default gro d.tif images are double-	oup and the group clicked in Explore	p to er Set File Associa	tion
Enter Multil Iser Mode	Configure	Hardware	Launch MDCStore	Loolo

26. Click OK

- 27. You may be asked to restart windows. Close all programs and save any documents before clicking **Yes**
- 28. The software has now been updated