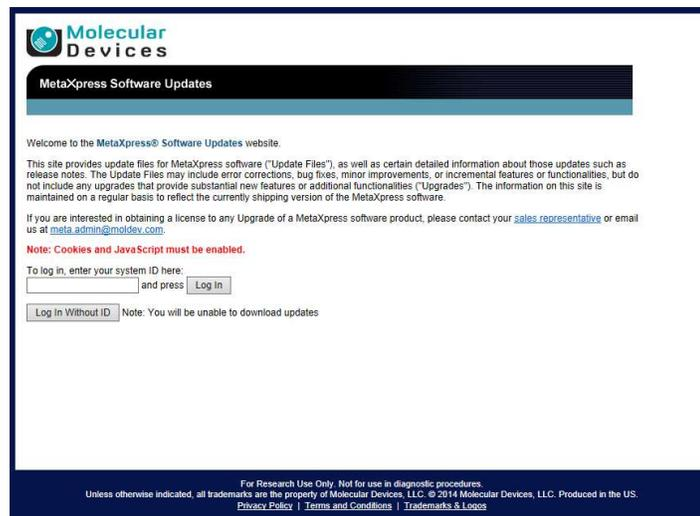


Updating MetaXpress® Software

1. Go to the MetaXpress Software Updates Page:
<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. In MetaXpress software on the main toolbar, click on File → About MetaXpress



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

Molecular Devices
MetaXpress Software Updates

[Main](#) **MetaXpress Software Updates**

Custom Module Editor

[License File Generator](#)

[2.5.13.3 Update](#)

[2.5 Release Notes](#)

[2.0.38.3 Update](#)

[2.0 Release Notes](#)

MetaXpress

[About MetaXpress](#)

[Code Generator](#)

[5.3.0.1 Update](#)

[5.3.0.0 Update](#)

[5.3 Release Notes](#)

Use the links at the left to view:

- **Release notes for MetaXpress Custom Module Editor 2.X update** 2 can be downloaded and installed, provided that prerequisites have been installed.
- **Release notes for MetaXpress 5.X updates.** Subsequent updates and applied, provided that a version of 5.X is already installed.
- **Release notes for MetaXpress 4.0 updates.** Subsequent updates applied, provided that a version of 4.0 is already installed.
- **Release notes for MetaXpress 3.1 updates.** Subsequent updates applied, provided that a version of 3.0 or 3.1 are already installed.
- **Release notes for MetaXpress 3.0 updates.** Subsequent updates applied, provided that a version of 3.0 is already installed.
- **Release notes for MetaXpress 2.0 updates.** Subsequent updates applied, provided that a version of 2.0 is already installed.
- **Release notes for MetaXpress 1.7 updates.** Subsequent updates applied, provided that a version of 1.7 is already installed.
- **Release notes for the latest MetaXpress 1.6 update.**

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 appropriate 32-bit or 64-bit 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

Molecular Devices
MetaXpress Software Updates

[Main](#) **MetaXpress 5.3.0.1 Update**

Custom Module Editor

[License File Generator](#)

[2.5.13.3 Update](#)

[2.5 Release Notes](#)

[2.0.38.3 Update](#)

[2.0 Release Notes](#)

MetaXpress

[About MetaXpress](#)

[Code Generator](#)

[5.3.0.1 Update](#)

[5.3.0.0 Update](#)

[5.3 Release Notes](#)

Posted 17 January, 2014

The MetaXpress 5.3.0.1 Update is available in both 32-bit and 64-bit versions.

The 32-bit version can be downloaded by clicking on the following link:

→ MetaXpress (150MB): [mx5_3_0_1_Win32_2014-01-15.zip](#)

If you have any previous 32-bit version of MetaXpress 5, then once you have the **Imaging Series Updater**. Point the Updater to the mx5_3_0_1_Win32_2014-01-15.zip

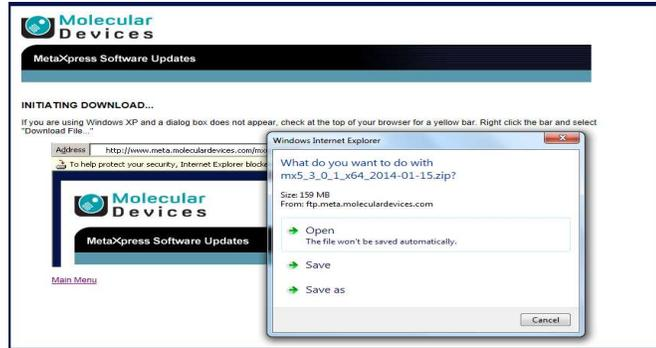
The 64-bit version can be downloaded by clicking on the following link:

→ MetaXpress (159MB): [mx5_3_0_1_x64_2014-01-15.zip](#)

If you have any previous 64-bit version of MetaXpress 5, then once you have the **Imaging Series Updater**. Point the Updater to the mx5_3_0_1_x64_2014-01-15.zip

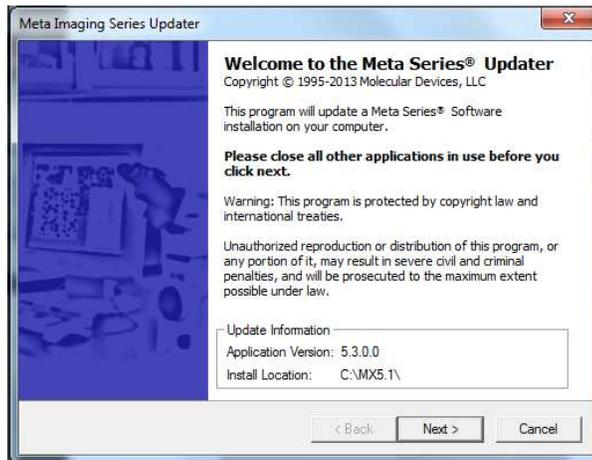
Note: The Updater communicates directly with our database to retrieve the latest update.

6. When the license agreement appears, click **I ACCEPT THE TERMS IN THIS AGREEMENT**
7. A bar or pop-up window will appear asking you if you want to open or save the file. Click on **Save As** and save the file to your hard drive in an easy to access location



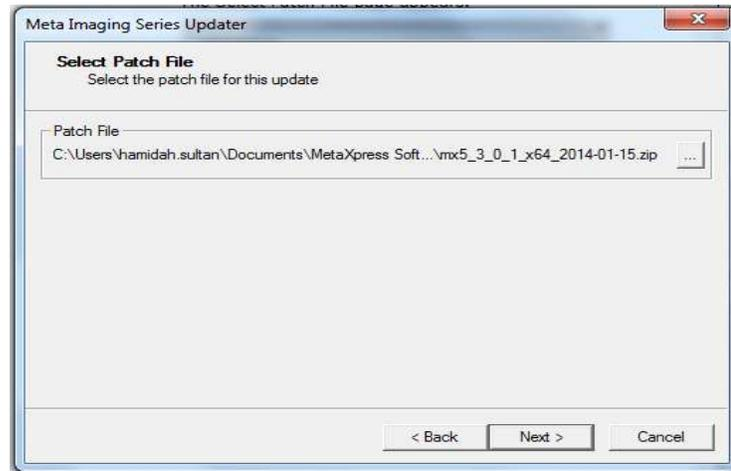
CAUTION! Do not attempt to unzip the update file. The MetaImaging Series® Updater unzips the file during the update process

8. From the Windows **Start** menu, select **Programs → MetaXpress 5.x → Meta Imaging Series Updater**



9. Click **Next**
10. Click on the ellipsis (...) button and the Windows Open dialog will appear
11. Locate the update file you just downloaded

12. Click **Next**



13. The **Hardware Key Confirmation** page will appear



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**

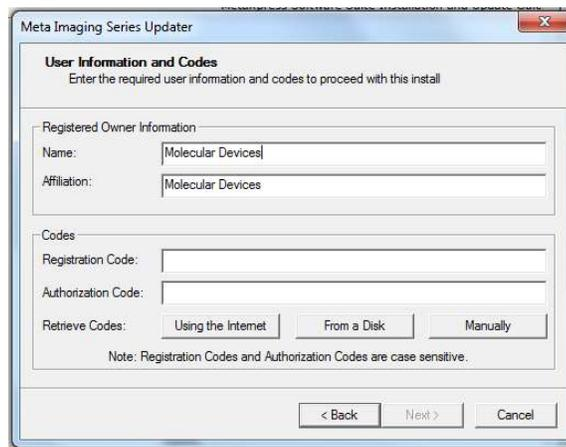
15. If you are using a network key to update the software on an analysis (offline) workstation computer, select **Network Key Server**, type the server address (static IP address or full computer name) where the network key is installed, and then click **Read System ID from Key**

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



The screenshot shows a Windows-style dialog box titled "Meta Imaging Series Updater". The main heading is "User Information and Codes" with the instruction "Enter the required user information and codes to proceed with this install". The dialog is divided into two sections: "Registered Owner Information" and "Codes". In the "Registered Owner Information" section, there are two text input fields: "Name:" containing "Molecular Devices" and "Affiliation:" containing "Molecular Devices". In the "Codes" section, there are two empty text input fields for "Registration Code:" and "Authorization Code:". Below these fields are three buttons labeled "Using the Internet", "From a Disk", and "Manually" under the heading "Retrieve Codes:". A note at the bottom of the dialog states "Note: Registration Codes and Authorization Codes are case sensitive." At the bottom of the dialog are three buttons: "< Back", "Next >", and "Cancel".

19. Enter your Name and Affiliation (optional)

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

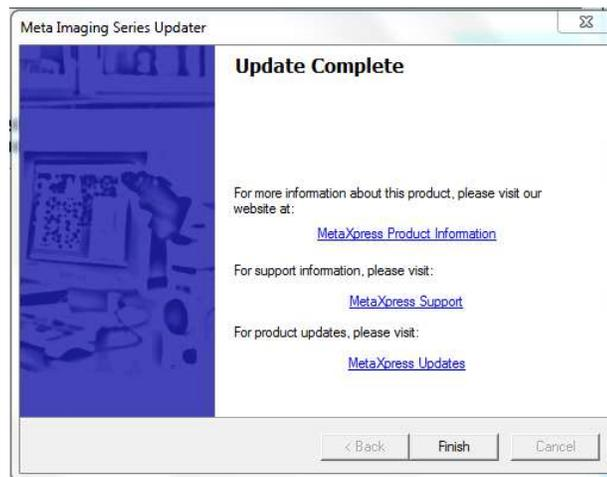
- a. If the computer is connected to the internet, click **Using the Internet**. The Updater retrieves the codes and displays them in the appropriate fields.
- b. If the codes are on a disk, insert the disk and click **From a Disk**. The Updater copies the codes from the disk to the appropriate fields.

- 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 <http://www.meta.moleculardevices.com/authorize/> . You will be asked to enter your System ID.

21. Click **Next**

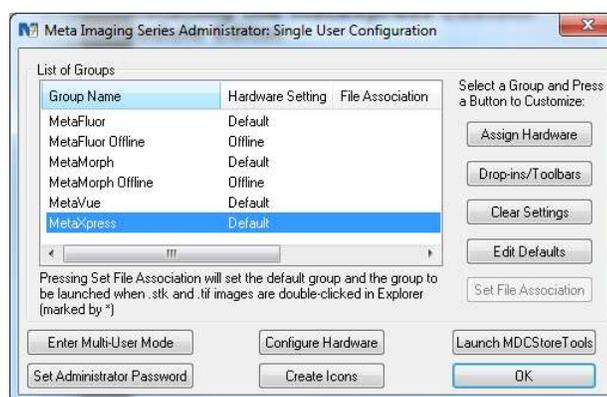
22. The **Ready to Proceed** dialog will appear. Click **Update**

23. When the installation is complete, the **Update Complete** screen appears



24. Click **Finish**

25. The Meta Imaging Series Updater opens the **Meta Imaging Series Administrator** screen



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