



MetaXpress[®] 6 FAQ

Drop-Ins and Custom Toolbars in MetaXpress

Date Revised 07/07/15 Version B



© 2012-2015. Trademarks property of Molecular Devices, LLC or their respective owners.
For research use only. Not for use in diagnostic procedures.

Frequently Asked Questions Overview

The purpose of this FAQ guide is to guide the user in enabling and disabling drop-ins as well as create custom toolbars for **MetaXpress** through the **Meta Imaging Series Administrator** software.

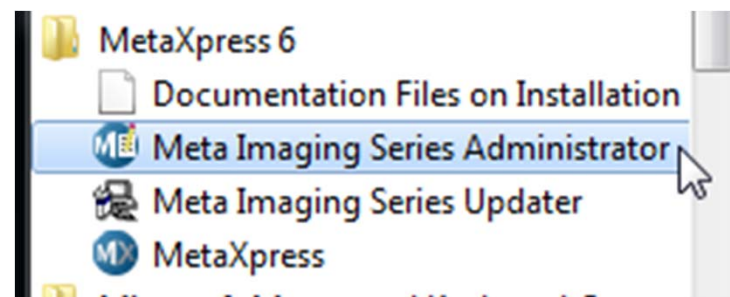


What is a Drop-in?

Many software functions in MetaXpress are enabled through drop-ins

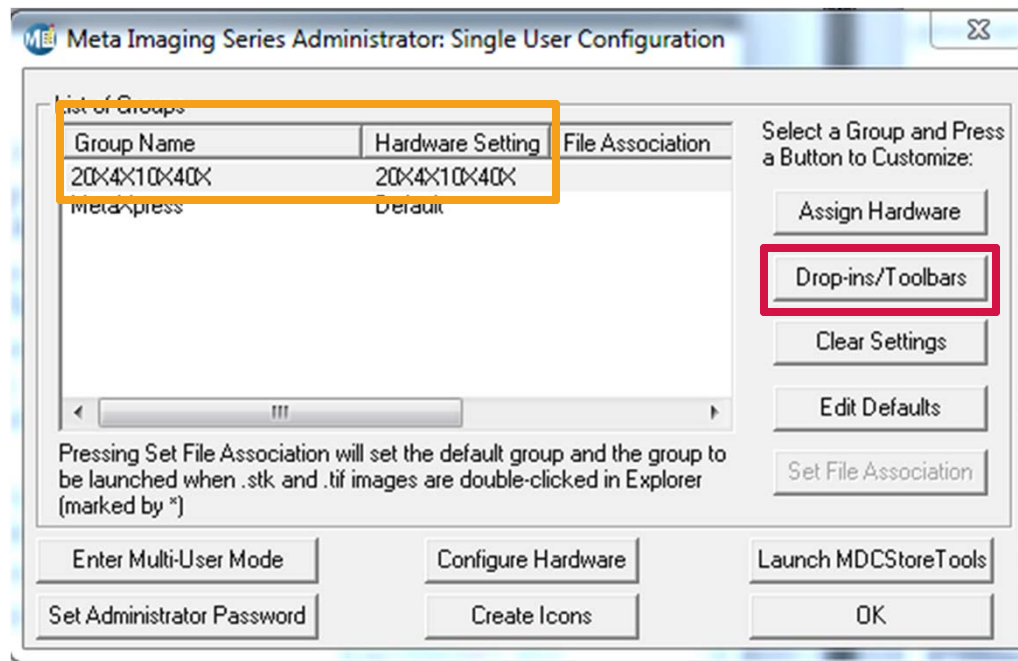
- Some drop-ins are standard and others need to be purchased (i.e. application modules, 3D Autoquant Deconvolution, etc.)
- The **Meta Imaging Series Administrator** software allows you to enable or disable these functions for each group (refer to corresponding chapters on creating groups and hardware profiles)

If you have just purchased an option, you must first update the software prior to enabling the Drop-in. Refer to corresponding chapter on updating MetaXpress.



Enable / Disable a Drop-in

1. Exit out of the **MetaXpress** software
2. Open the **Meta Imaging Series Administrator** software
3. Highlight the group in which you would like to enable or disable the Drop-in
4. Click on the **Drop-ins/Toolbars** button

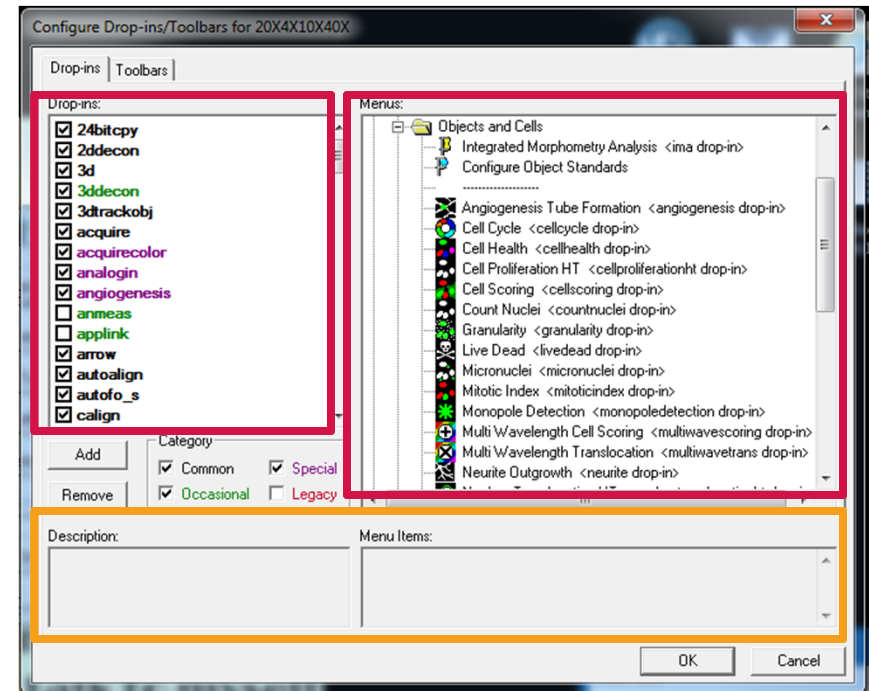


Enable / Disable a Drop-in

5. Select the **Drop-ins** tab
6. Select or deselect the check box next to the appropriate Drop-in in the left side panel
7. If you are unsure of the function, highlight the Drop-in and a description will appear at the bottom of the dialog

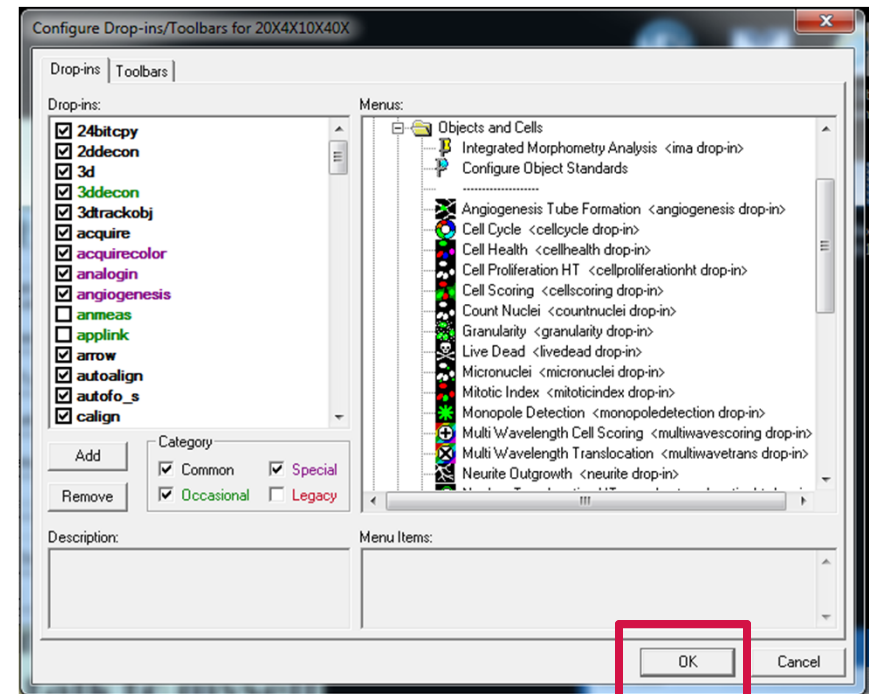
NOTE If the drop-in is not visible, enable all check boxes in the **Categories** section

8. An enabled Drop-in will appear on the right side panel under the main menu category found in MetaXpress



Enable / Disable a Drop-in

10. Click **OK** to close the **Configure Drop-ins/Toolbars** dialog
11. Repeat steps 3-8 to enable or disable the Drop-in for other groups
12. Click **OK** to exit the **Meta Imaging Series Administrator** software
13. Open MetaXpress to verify the Drop-in has been enabled/disabled



List of Common Drop-ins

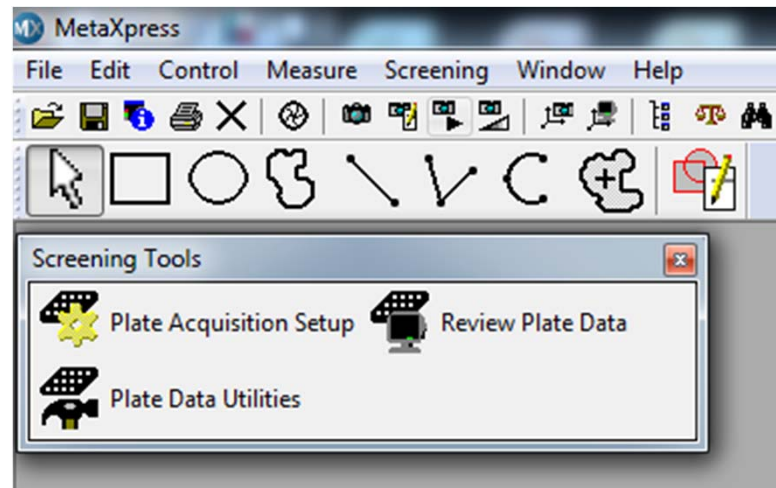
Function name	Drop-in name in Meta Imaging Series Administrator
Angiogenesis Tube formation	angiogenesis
Cell Cycle	cellcycle
Cell Health	cellhealth
Cell Proliferation HT	cellproliferationht
Cell Scoring	cellscoring
Count Nuclei	countnuclei
Granularity	granularity
Live Dead	livedead
MicroNuclei	micronuclei
Mitotic Index	mitoticindex
Monopole Detection	monopoledetection
Multi Wavelength Cell Scoring	multiwavescoring
Multi Wavelength Translocation	multiwavetrans
Neurite Outgrowth	neurite
Nuclear Translocation HT	nucleartranslocationht
Transfluor	gpcr
Transfluor HT	transfluorht
Translocation	translocation
Translocation Enhanced	translocationenhanced
MDMA	3dtrackobj



What is a Toolbar?

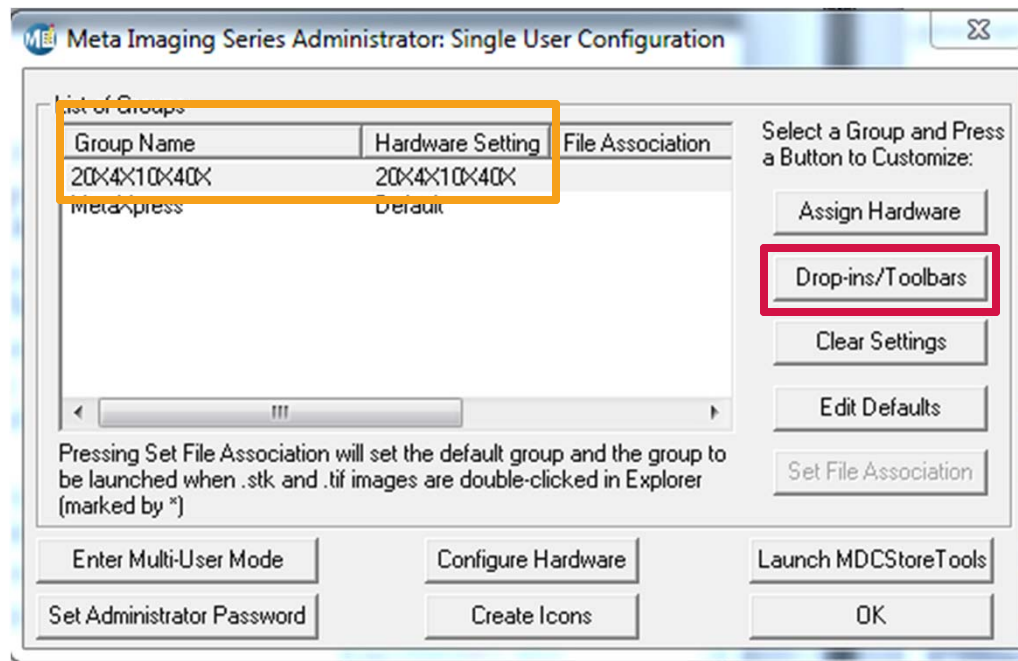
Toolbars in MetaXpress are a collection of menu items, functions, and journals that can be configured to appear under the main menu or as floating boxes

- Toolbars can be customized with user defined text and icons
- The **Meta Imaging Series Administrator** software allows you create toolbars for each group (refer to corresponding chapters on creating groups and hardware profiles)



Creating a Toolbar

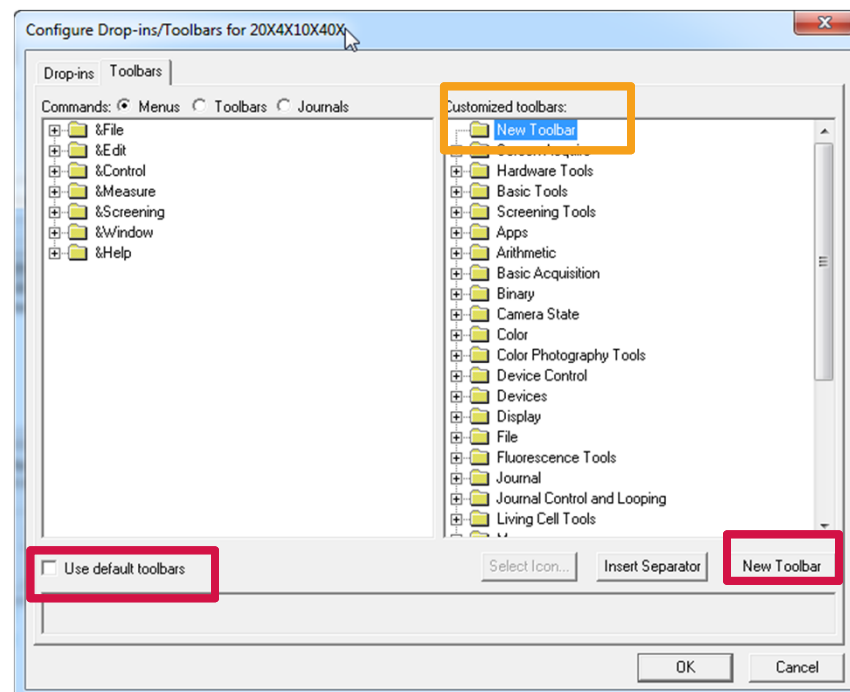
1. Exit out of the **MetaXpress** software
2. Open the **Meta Imaging Series Administrator** software
3. Highlight the group in which you would like to create a toolbar
4. Click on the **Drop-ins/Toolbars** button



Creating a Toolbar

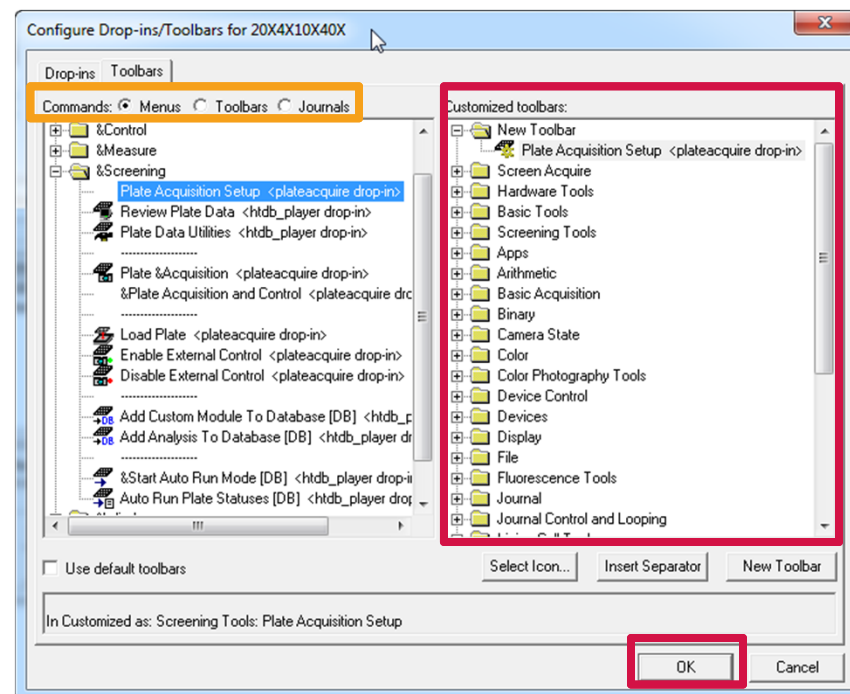
5. Select the **Toolbars** tab
6. Uncheck the **Use default toolbars** check box
7. Click on the **New Toolbar** button
8. This will create a new folder called “New Toolbar” in the right side panel
9. Right-click and select **Rename** or double click on the folder to edit the name

NOTE Remember this name as you will need it when enabling the toolbar in MetaXpress



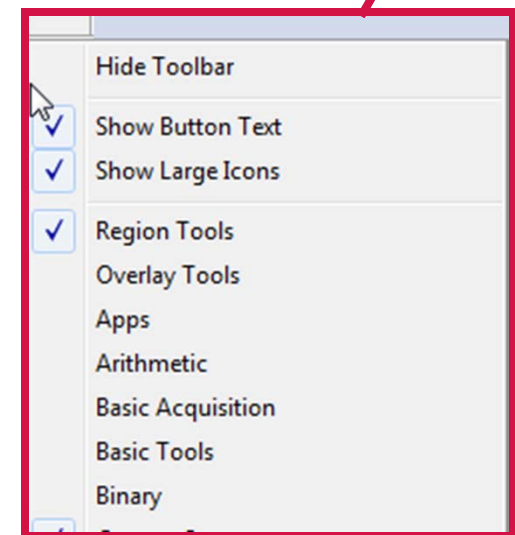
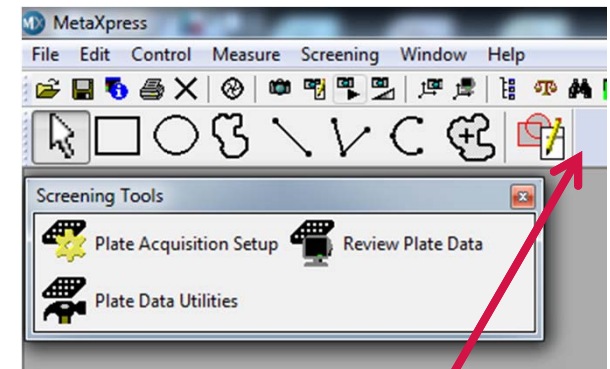
Creating a Toolbar

- Expand the folders in the left side panel to display menu items, toolbars, and journals
- Drag and drop items from the left side to the right side panel under the folder that was just created
- Right-click on the item to edit the text that will appear in MetaXpress
- Click on the **Select Icon** button to assign a different icon
- Change the selections next to **Commands** for more options



Creating a Toolbar

15. When you have finished adding and configuring items, click on the **OK** button
16. Repeat steps 3-15 for additional groups
17. Click **OK** to exit the **Meta Imaging Series Administrator** software
18. Open MetaXpress
19. Right click anywhere beneath the main menu and enable the new toolbar you just created
20. Right click on the toolbar itself to enable display button text and change icon size
21. Drag and drop the toolbar to configure its location or to create a floating window



Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- User Forum: <http://metamorph.moleculardevices.com/forum/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments





ADVANCING PROTEIN AND CELL BIOLOGY