



ImageXpress Taskbar Release Notes

Version 6.7

New Features in IX Taskbar 6.7

- Support added for Confocal HT.ai instrument.
 - Updated shading plate and GP slide selections
 - Updated following workflows to reduce likelihood of damaging the target:
 - Verify A1 Center
 - Measure Pixel Sizes
 - Parfocality and XY Offsets
 - Set up Shading Correction
 - Adjust Stage Position journal: Added option to adjust intensity level
 - New Reset Interlocks button added to Run a Plate, Scan Slides, and System Maintenance taskbars.

Improvements and Bug Fixes 6.7.1

- IX Shutdown Journal: No longer prompts twice to open the door.
- Set up Shading Correction: Improved workflow
 - Allows user to save widefield images with confocal naming
 - Allows user to specify combinations in advance for slide shading correction
- Improved Verify A1 Center workflow
 - Allows user to adjust the focus and exposure time
- Updated simplified menu

New Features in IX Taskbar 6.6

- Support added for Water Immersion objectives
- Updated custom menu with new Water Immersion dialogs
- In Measure Pixel Sizes, added new option (test montage) for measuring/adjusting pixel sizes.

Improvements and Bug Fixes 6.6.1

- Updated Set up Slide Dimensions to prevent light exposure while performing visual measurement inside IXM; close door automatically after performing visual measurements.
- Shading correction log now appends always instead of offering overwrite option.
- Improved messaging for Adjust Correction Collar
- Updated IX Shutdown Journal: If WI objective is last objective selected, always open door and prompt user to remove plate
- Updated Close Door and Load Plate Journals to clear device errors (for laser safety interlocked systems).
- Updated Set up Shading Correction: Air objectives are done before WI objectives. Slide shading correction now has correct confocal information in image names and removes special characters from the image names.
- Updated Parfocality and XY Offsets, Measure Pixel Sizes to prompt user whether to open door for GP-2 slide, corrected color plastic slide labels in user prompt. Run all air objectives prior to water objectives.
- Updated Slide Autofocus Wizard: Prompt user whether to open door or not. Run all air objectives prior to water objectives.
- Updated Laser Autofocus on Slide to add compatibility for Water Immersion objectives, added more retries, limited single search ranges for SWD objectives to reduce chance of collision.
- Minor bug fixes in Create Scan Areas

New Features in IX Taskbar 6.2

- Support added for ImageXpress Nano system with MetaXpress software.
- Set up Shading Correction now supports creating shading images in confocal as well as widefield modes on an ImageXpress Micro Confocal system.
- Taskbar now works if installed in MetaXpress installation folder that is not a top-level folder on the drive.
- Slide scanning tools now compatible with custom slide holders.
- New lookup table makes it simpler to customize slide laser autofocus settings for the taskbar.
- Close Log button added to Analyze Images taskbar.
- Scale Images to Full Range and AutoScale Images now apply to all open images.
- Load Stack – Current Site added to Analyze Images taskbar.
- Load Montage – Current Well added to Analyze Images taskbar.
- Copy Mask and Add to Stack added to Analyze Images taskbar.
- IX Shutdown journal provided to turn off Transmitted Lamp (if system has Transmitted Light option) and prompt user to remove plate from system.
- Turn off Transmitted Lamp button added to Run a Plate taskbar if system has Transmitted Light option.
- Alternative routine for Parfocality only added to Parfocality and XY Offsets function.
- IX Startup journal has additional option to check liquid levels during startup and display warning if low (if system has Environmental Control option).

Improvements and Bug Fixes 6.2.3

- Laser autofocus on slide and related journals no longer give laser intensity error with MetaXpress 6.5.3 or 2x objectives.
- Create Scan Areas no longer shows MetaMorph cannot load images of this format error.
- Create Scan Areas now only offers custom FOV option if custom FOV is < 90%.
- A warning is now displayed if Live mode is on when Open Door or Eject Plate is selected.
- The optional objective verification step in the Startup journal now labels the "Closest" and "Farthest" objectives correctly.

Improvements and Bug Fixes 6.2.2

- Taskbar installer now correctly prompts installation of Center on Click function if not enabled already.

Improvements and Bug Fixes 6.2.1

- Custom menu updated.
- Eject Filter Cubes now moves the filter cube changer to the eject position as expected for widefield systems.
- Create Scan Areas now allows you to update an existing plate file.
- Corrected camera settings for ImageXpress Micro Confocal system.
- Simplified entry of names when using Perform Preview Scan and Create Scan Areas.
- Perform Preview Scan always prompts user to adjust/confirm exposure time and focus.
- Perform Preview Scan uses the correct wavelength selection chosen by user during setup.
- Set up Shading Correction for Slides allows you to run multiple filter/objective combinations without error.
- Set up Slide Dimensions improved when using visual measurement of slide in system.
- Improved default starting XY positions when calibrating A1 Center (running Verify A1 Center in Maintenance Mode).
- Create Scan Areas offers user option to retain Z series settings when updating an existing protocol.
- Create Scan Areas displays an error message if region selection requires more sites than available in Plate Acquisition Setup.
- Create Scan Areas provides more information to user when choosing between full field of view and custom field of view.
- Taskbar warns user that toolbar icons may be missing if installing custom menu and MetaXpress is not in default directory C:\MX6
- Parfocality and XY Offsets now provides the hardware settings name in the instructions provided to user for entering new parfocality values.
- Set up Shading Correction now drops the objective to a low Z position when changing objectives.
- Taskbar buttons rearranged for more logical workflow.
- Laser Autofocus on Slide function now shows an informative error message if an oil-immersion objective is selected.
- Create Scan Areas adds an "After Each Image" journal for the Plate Acquisition protocol to ensure that acquired images have an appropriate stage label assigned to them.
- Image showing fluorescent plastic GP slides has updated labeling matching the slides in the accessory kit.