



ImageXpress Taskbar Release Notes

Version 6.2

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.