



MetaXpress[®] 6 Software Guide

IXM Taskbar Installation and Overview

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.

Chapter Purpose

The purpose of this chapter is to guide the user in installing the **IXM Taskbar** and give a brief overview.

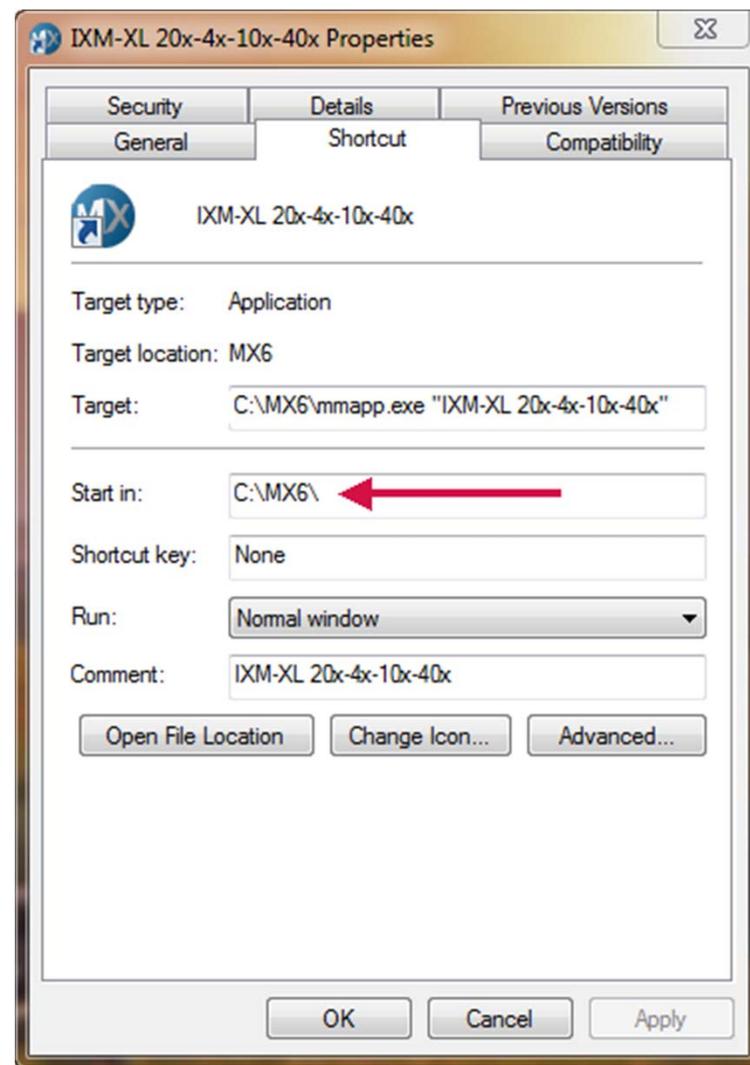
The **IXM Taskbar** is a collection of tools intended to enhance and streamline common imaging workflows. For details on specific taskbar workflows, refer to corresponding chapters and FAQ.

The **IXM Taskbar** is available on the User Forum:
<http://metamorph.moleculardevices.com/forum/>



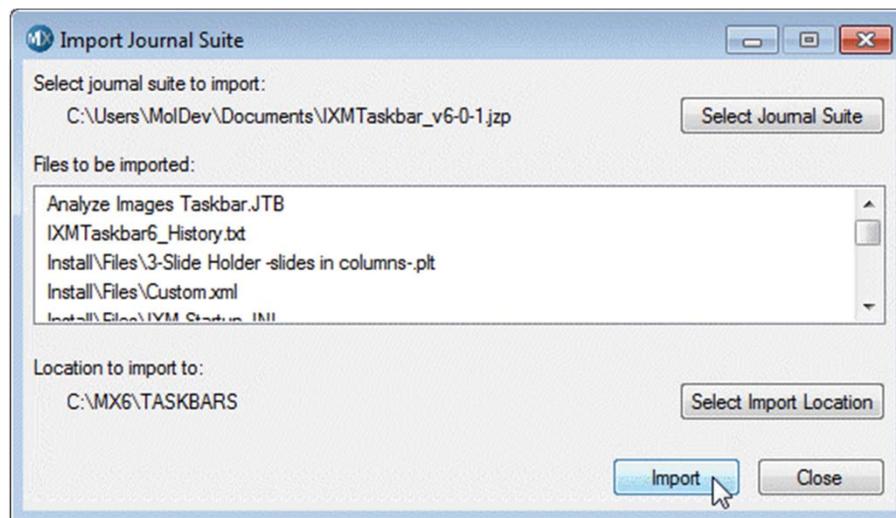
Installing the IXM Taskbar: Verify MX Folder

1. Identify the main folder where MetaXpress 6 is installed.
 - Default folder location is usually C:\MX6\
2. To verify the MX6 folder
 - Right-click on any of the MetaXpress shortcut icons and select Properties
 - Note the location listed in **Start in:**
3. Open the folder identified in Step 2 and create a new folder called **Taskbars** if it does not exist already
 - Molecular Devices recommends to delete all files in the **\Taskbars** and **\Startup** folders
4. Power on the system and start the MetaXpress software (any hardware profile)



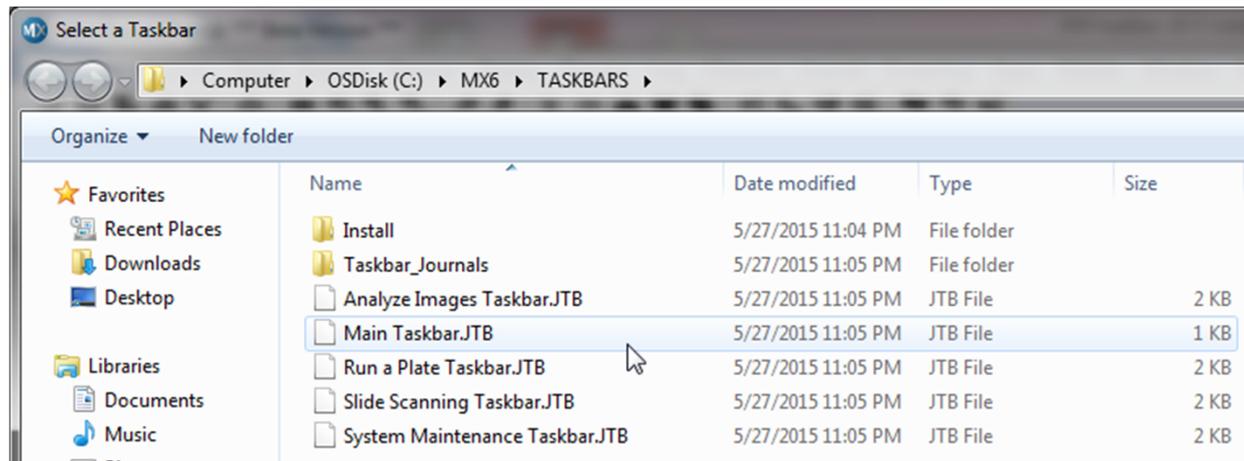
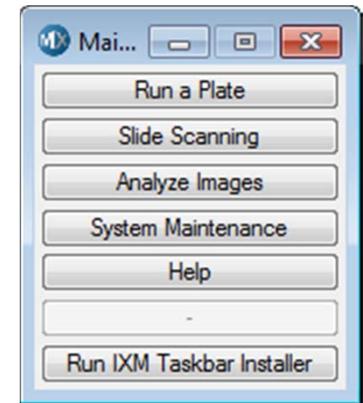
Installing the IXM Taskbar: Import Journal Suite

- Download the latest version of the IXM Taskbar from the User Forum (<http://metamorph.moleculardevices.com/forum/>). If the file is a .zip file, unzip the folder to for the .jzp file needed in step 6 below.
- Import the journal suite (.jzp):
 - In the main menu, select **Control > Journal > Import Journal Suite**
- Click on **Select Journal Suite** and locate the .jzp file
- Click on **Select Import Location** and browse to the **\Taskbars** subfolder in the MX6 folder
 - This is usually **C:\MX6\Taskbars**
- Click **Import**
- Click on Close when complete

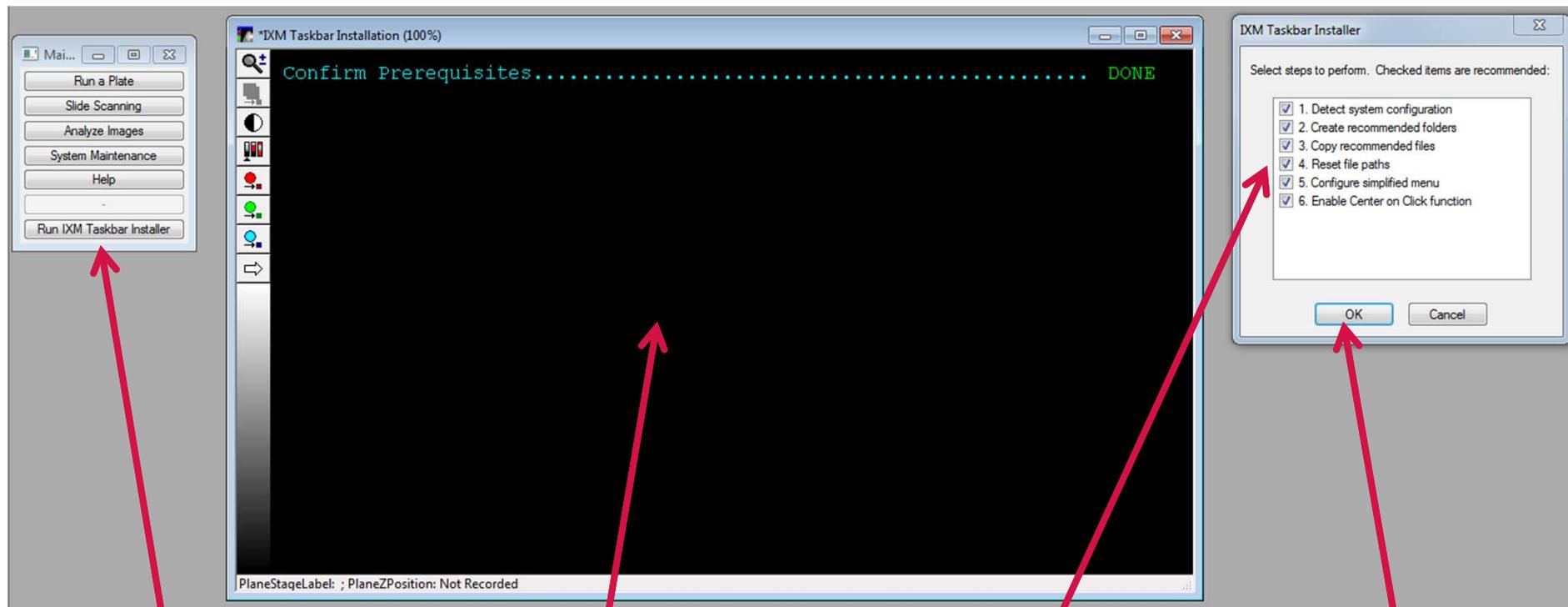


Installing the IXM Taskbar: Load Taskbar

11. In the main menu, select **Control > Journal > Taskbars > Load Taskbar**
12. Select the **Main Taskbar.JTB** file in the \Taskbar subfolder of the MX6 folder
 - This is usually **C:\MX6\Taskbars**
13. The **Main Taskbar** should now appear in MetaXpress



Installing the IXM Taskbar: Run Taskbar Installer



14. Click on **Run IXM Taskbar Installer**

15. Status window appears

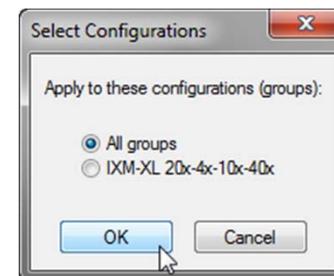
16. Select steps to perform

17. Click **OK**



Installing the IXM Taskbar: Run Taskbar Installer

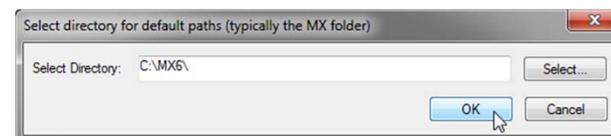
18. If prompted, select current group (configuration) or **All Groups** to install and apply configurations. Molecular Devices suggests to install for all groups. Click **OK**.



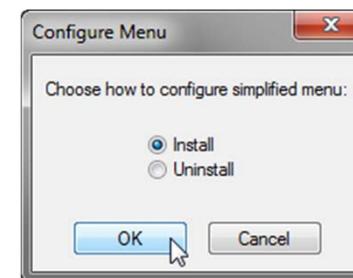
19. If prompted, select current user or **All users** to install and apply configurations. Click **OK**.



20. When prompted for default paths, choose the base folder for MetaXpress (usually **C:\MX6**). Click **OK**.

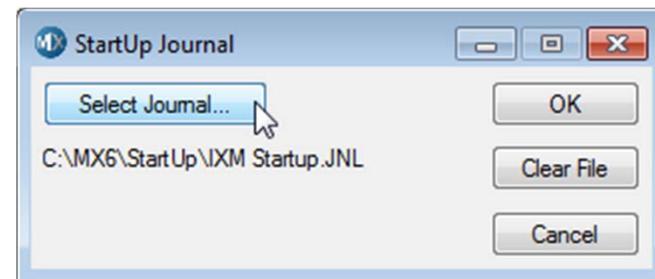


21. Choose to install or uninstall the simplified menu. Refer to corresponding chapters on simple vs. default menu for more details.



Installing the IXM Taskbar: Startup Journal

22. (OPTIONAL) MetaXpress can be configured to automatically display the IXM Taskbar upon restarting if it is accidentally closed
23. In the main menu, select **Control > Journal Control > Startup Journal**
24. Click on **Select Journal**
25. Browse for the IXM Startup.JNL file located in the \Startup subfolder in the MX6 folder (usually **C:\MX6\Startup**)
26. Click **OK**
27. Steps 22-26 will need to be repeated for each hardware profile.



Installing the IXM Taskbar: Restart MetaXpress

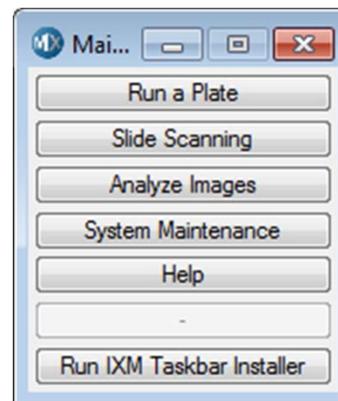
```
*IXM Taskbar Installation (100%)  
Confirm Prerequisites..... DONE  
1. Detect system configuration..... DONE  
2. Create recommended folders..... DONE  
3. Copy recommended files..... DONE  
4. Reset file paths..... DONE  
5. Install simplified menu..... DONE  
6. Enable Center on Click function..... DONE  
  
Recommended:  
Go to Journal > Journal Control > Startup Journal  
Select C:\MX6\StartUp\IXM Startup.JNL and click OK  
  
To complete the install, exit and restart MetaXpress  
  
PlaneStageLabel: ; PlaneZPosition: Not Recorded
```

28. If prompted, restart MetaXpress



Taskbar Overview: Main Taskbar

- Links to additional taskbars
 - **Run a Plate**
 - **Slide Scanning**
 - **Analyze Images**
 - **System Maintenance**
- Help (on all taskbars)
 - How to obtain support
 - Taskbar Version
 - System ID
 - Group (configuration/profile)



Taskbar Overview: Run a Plate Taskbar

Open and Close Door to access plates

Links to common plate tasks

- Plate Acquisition Setup
- Load/Save Protocol
- Acquire Plate

Adjust Correction Collar

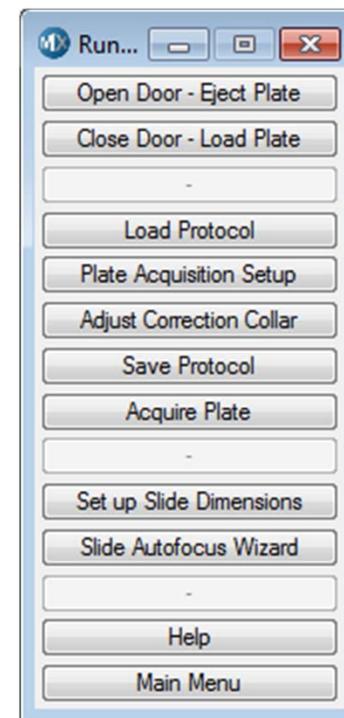
- Access objective to adjust collar for current plate type
- Usually found on 20x and higher magnification

Set up Slide Dimensions

- Recommended for slides with regularly spaced features
- Guided set up of a plate file for a slide

Slide Autofocus Wizard

- Create autofocus settings for slides



Taskbar Overview: Slide Scanning Taskbar

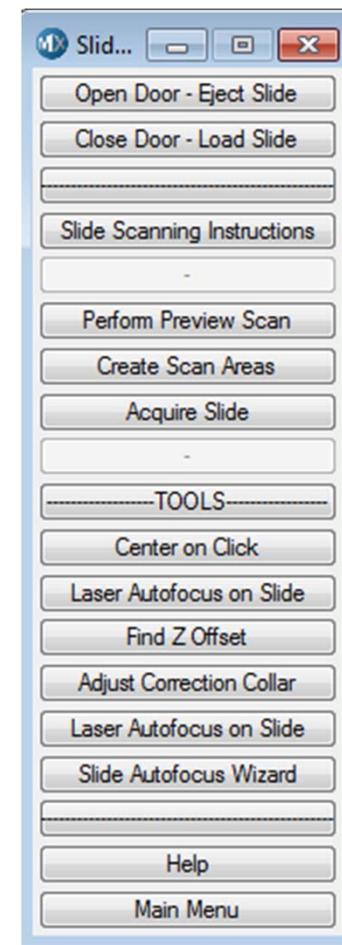
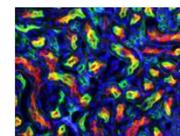
Open and Close Door to access slides

Slide Scanning Workflow for slides with irregularly spaced features (e.g. tissue slice/TMA)

- **Perform Preview Scan** (run low-magnification scan)
- **Create Scan Areas** (define regions of interest)
- **Acquire Slide** (acquire regions at high magnification)

Slide Scanning tools to use during setup

- **Center on Click**: move to a region of an image by clicking on it
- **Slide Autofocus Wizard**: create autofocus settings for slides
- **Adjust Correction Collar**: access objective to adjust collar for current slide



Taskbar Overview: Analyze Images Taskbar

Select Current Plate

- Selects the last acquired plate in Review Plate Data

Image Scaling (display setting)

- Scale Image to Full Range
- **AutoScale Image**

Measure Images to Estimate Module Settings

- **Estimate Module Settings**: quick tool to calculate width, area, intensity
- **Calipers**: measure width
- **Linescan**: view object and background intensities

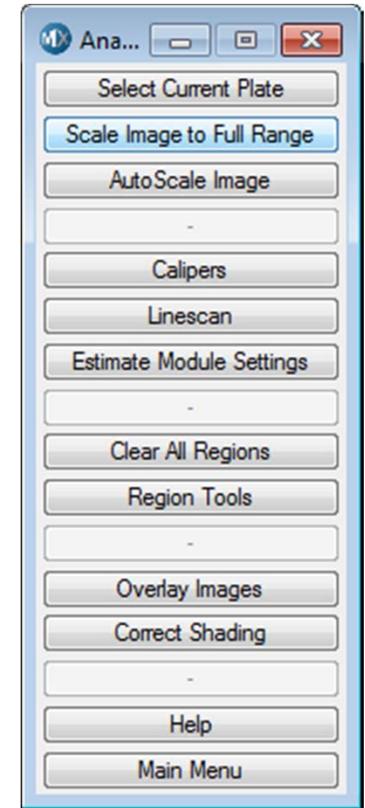
Regions

- **Clear All Regions**
- **Region Tools**: Opens the Region Tools toolbar (if not already open)

Overlay Images

Correct Shading

- Apply shading correction to an already acquired image



Taskbar Overview: System Maintenance Taskbar

Open and Close Door

- Option to leave the sample stage in place

Access objectives and filters

- **Adjust Correction Collar**
- **Eject Filter Cubes**

System calibration

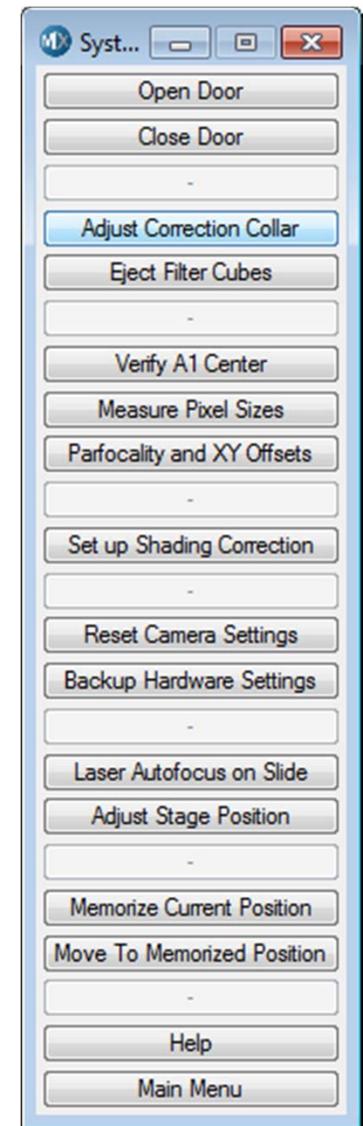
- **Verify A1 Center**: check the XY calibration using the slide holder
- **Measure Pixel Sizes**: verify the objective calibrations
- **Parfocality and XY Offsets**: measure objective offsets

Set up Shading Correction

- Automatically create shading correction images for plates or slides

Additional Tools

- **Reset Camera Settings** to factory defaults
- **Backup Hardware Settings** (illumination, magnification, calibration)
- **Laser Autofocus** on a slide
- Interactively **Adjust Stage Position**
- Memorize and **Move to a Stage Position**



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