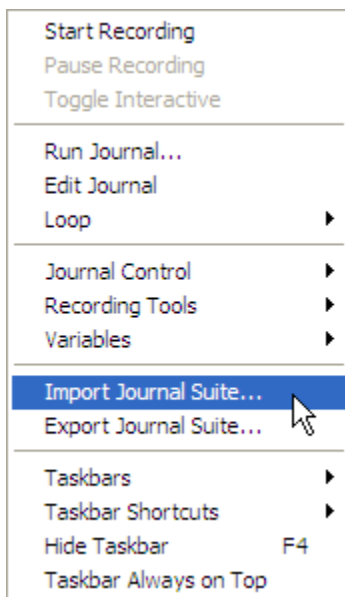
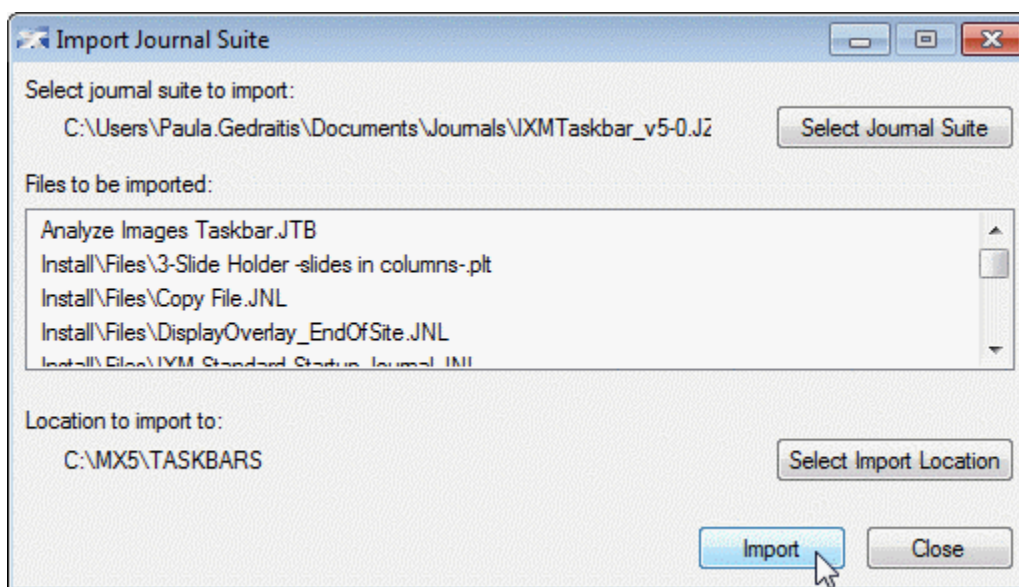


IXM Taskbar version 5.0 - Installation

1. Copy the provided IXMTaskbar_v5-0.jzp file to the IXM computer. The IXM must be using MX 4.0 or MX5.0.
2. Start up MetaXpress and go to **Journal > Import Journal Suite**.

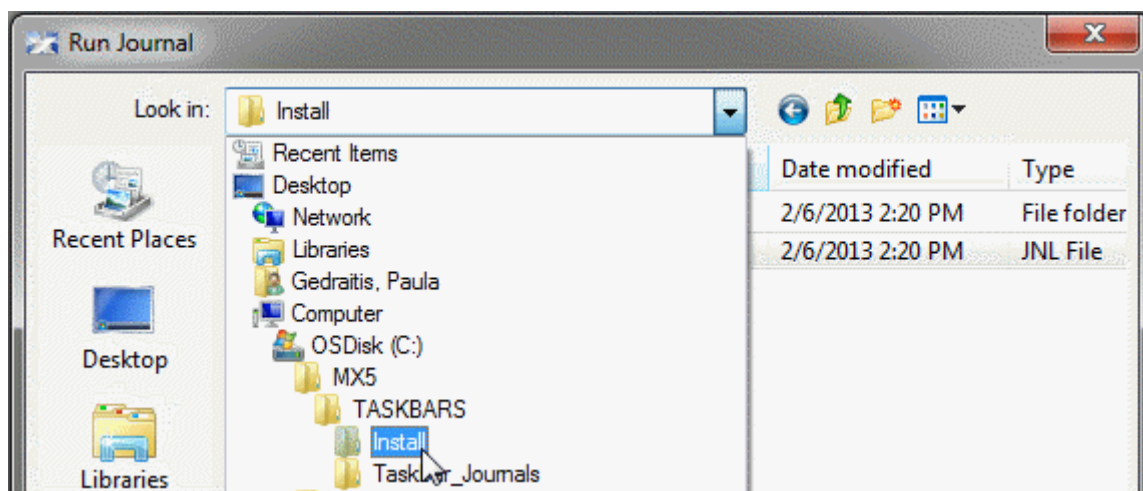
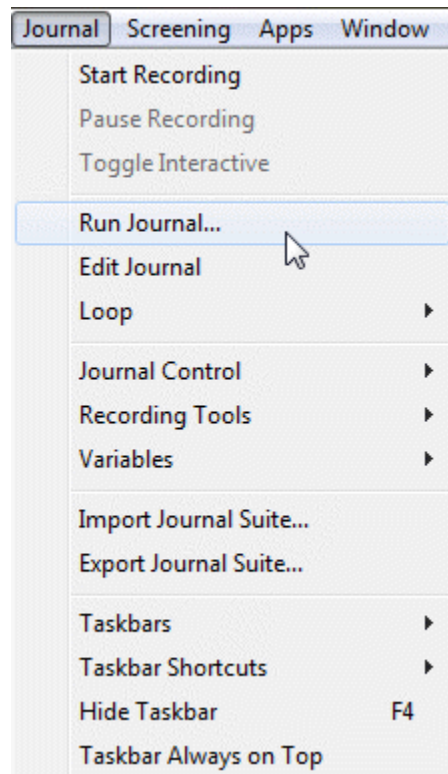


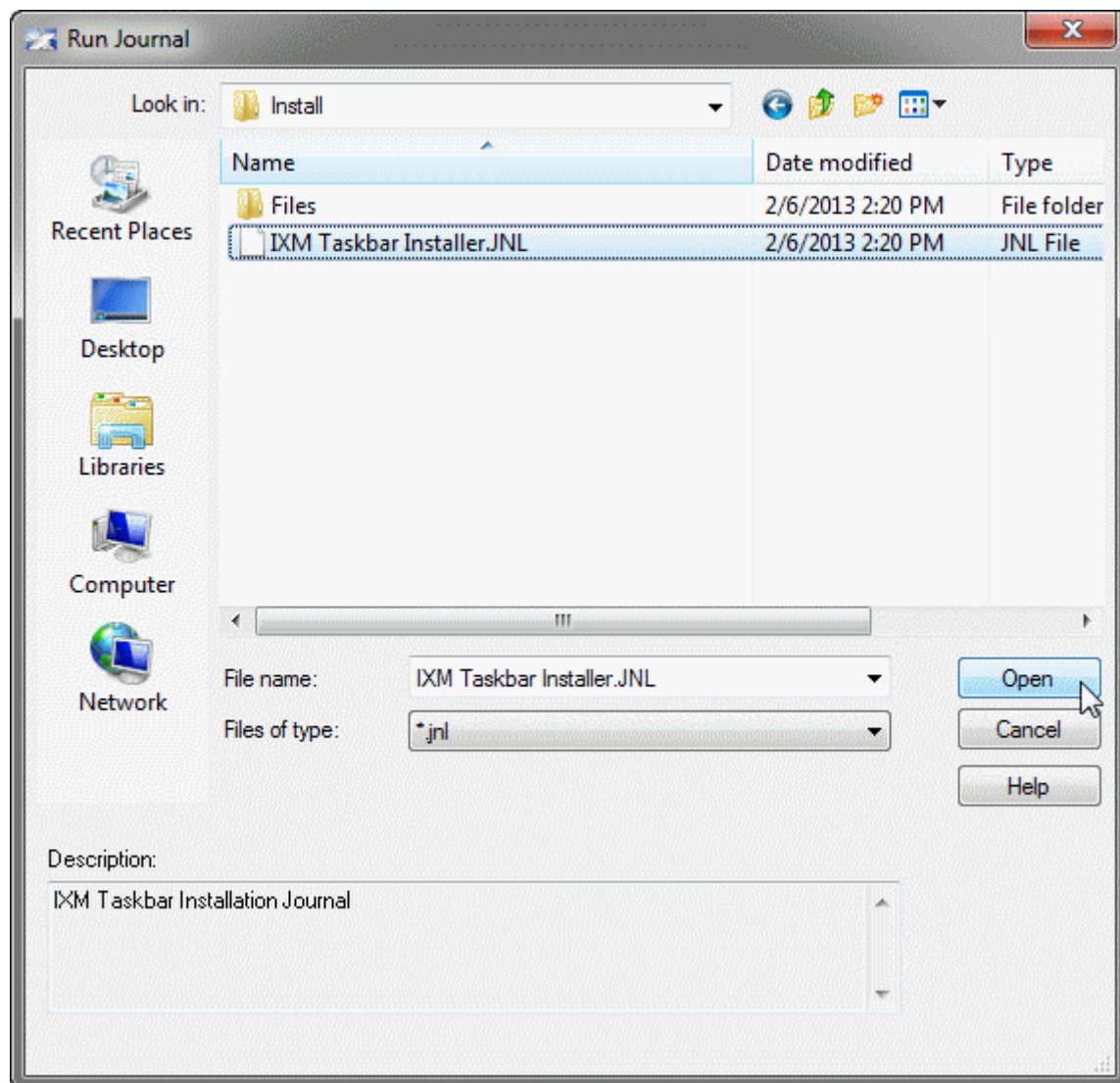
3. Select the file you copied over, and the **C:\MX4\Taskbars** directory or **C:\MX5\Taskbars** directory (if this directory does not exist in the current MetaXpress folder, please create it). Click **Import**.



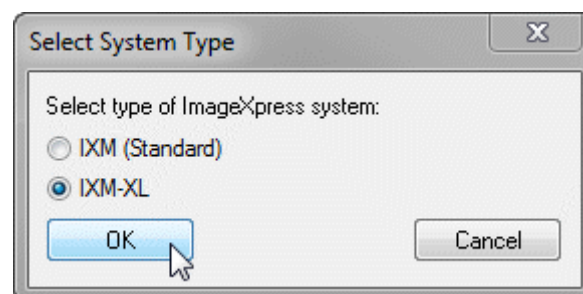
NOTE: The following steps must be repeated for every group (configuration).

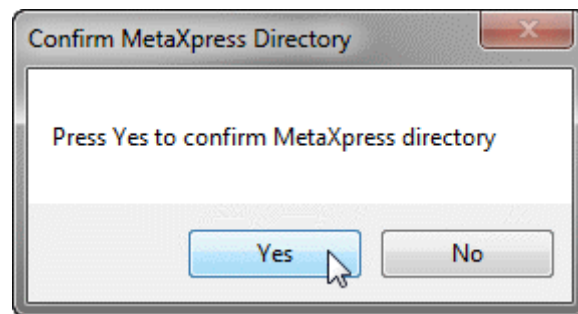
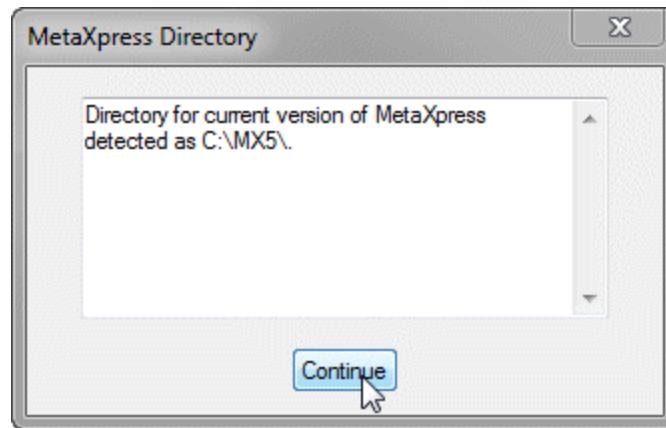
4. Go to **Journal > Run Journal**, navigate to the **C:\MX5\Taskbars\Install** folder and select the **IXM Taskbar Installer** journal.



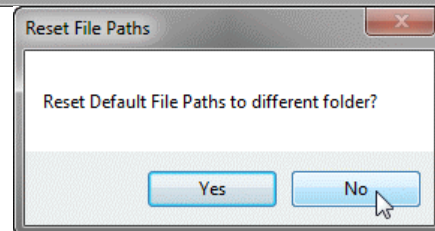
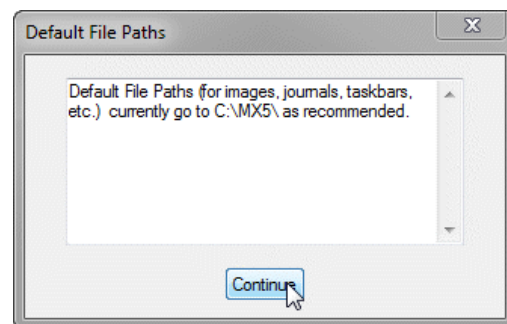
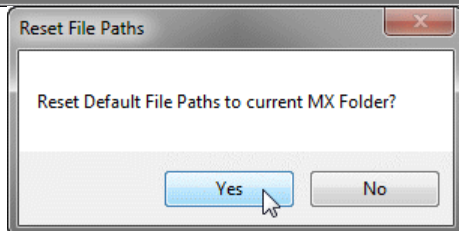
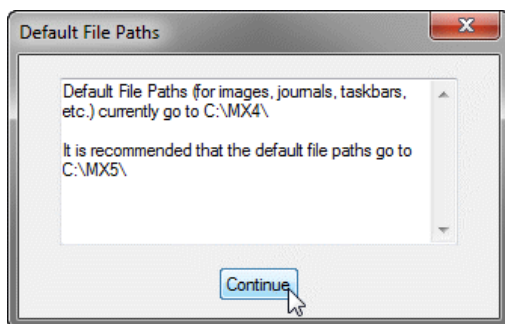


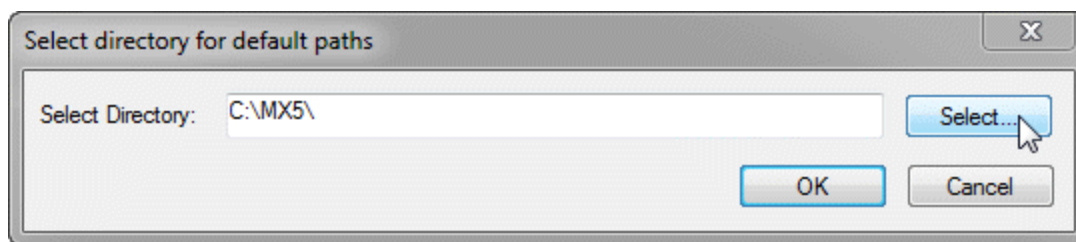
5. Follow the prompts to select the type of IXM (Standard or XL) and to confirm the directory that the current version of MX is installed in (typically C:\MX4\ or C:\MX5\).



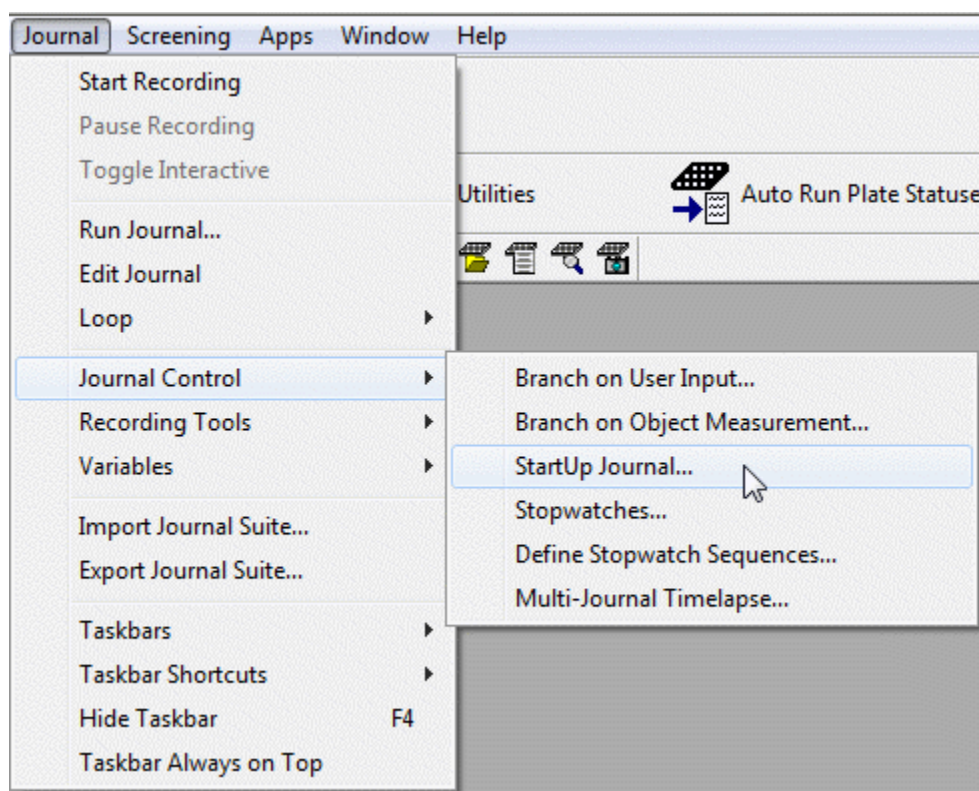
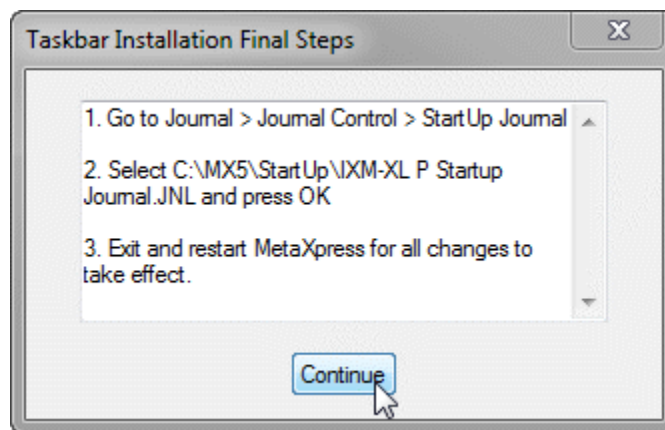


6. If the default file paths (for journals, images, data, etc.) do not point to the current MX folder, you will be provided with an option to reset them to the current folder.

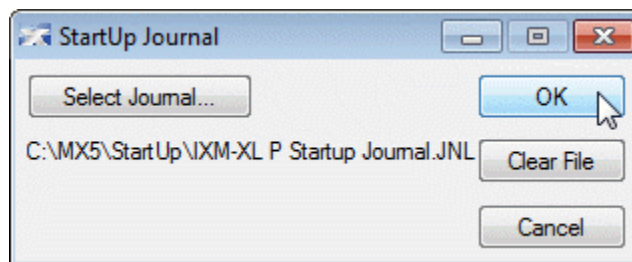
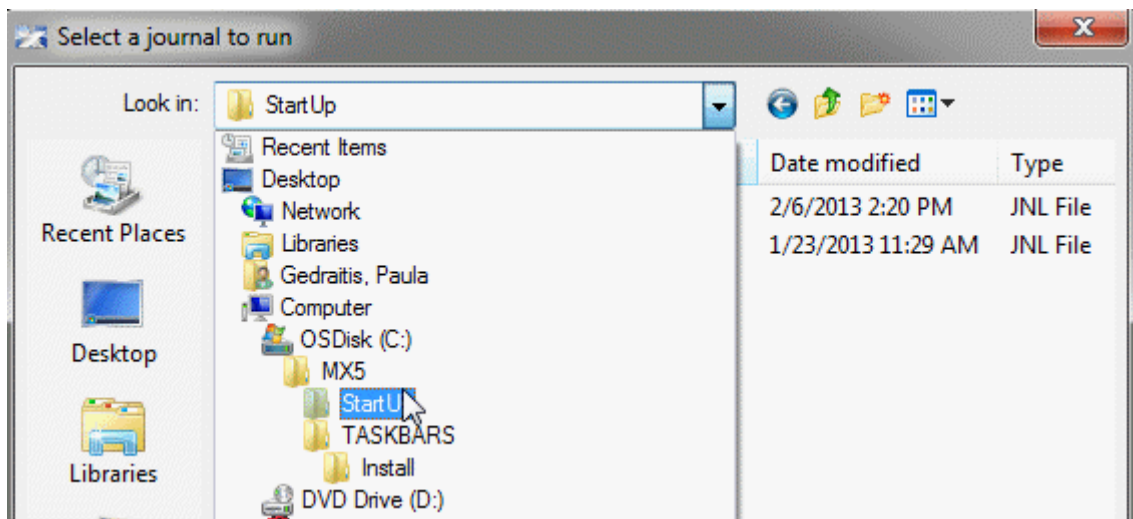
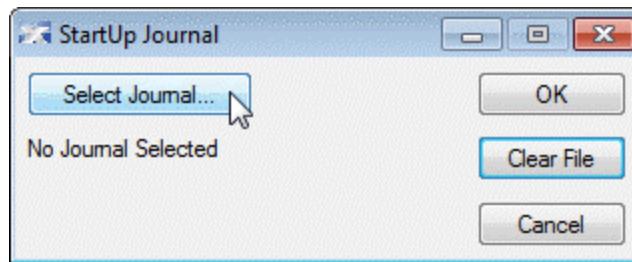




7. When the install journal has completed, you will be prompted to set the startup journal (see below). Go to **Journal > Journal Control > StartUp Journal**.

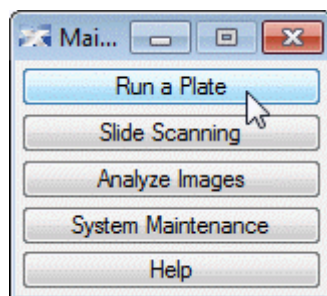


8. Press **Select Journal** and navigate to the **C:\MX4\StartUp** directory or **C:\MX5\StartUp** directory. Select the startup journal found there. It should be the appropriate journal for the IXM system, either Standard or XL. Press OK.

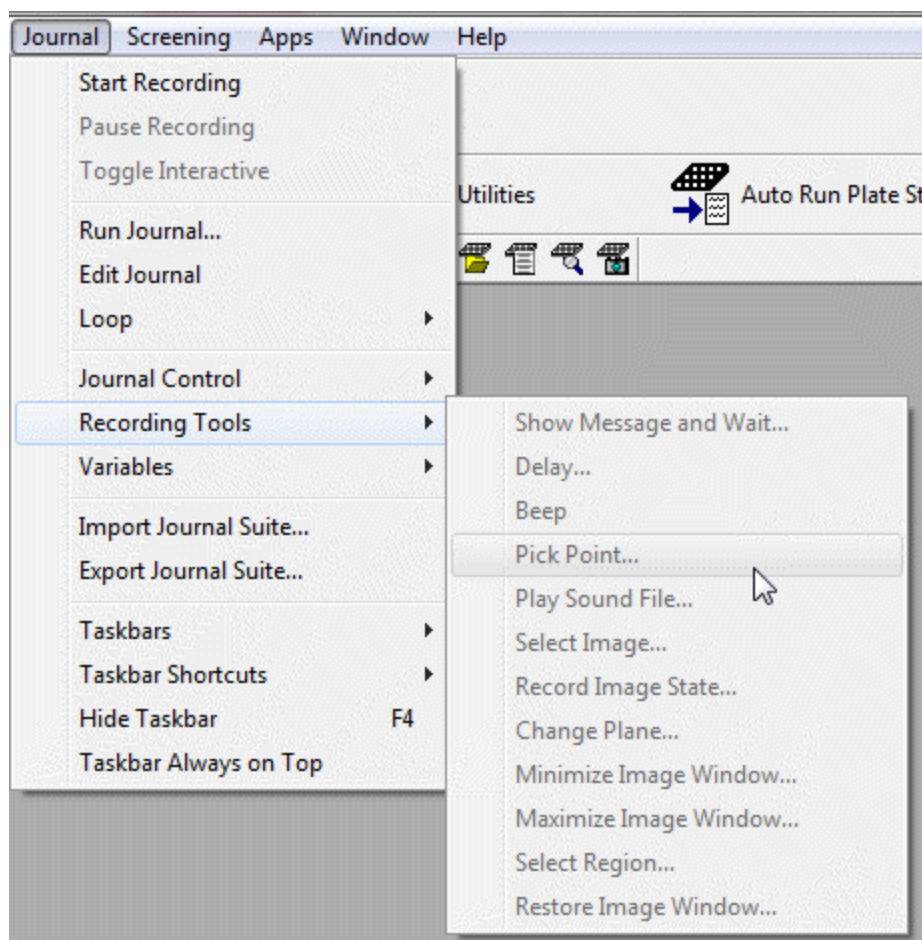


9. Now exit and restart MetaXpress for changes to take effect.

10. Verify that the taskbar is loading properly.



11. Also verify that Pick Point appears in the Journal > Recording Tools menu (it may be grayed out, which is fine).



12. Repeat steps 4-11 for all groups (configurations).

NOTE: This installation is easier if you perform it prior to creating additional groups (configurations) in Meta Imaging Series Administrator. Additional groups should then be created using the option to Copy Settings from an existing group.

Please contact Molecular Devices Tech Support with any issues or questions.