

MetaXpress Analysis Taskbar Installation Guide

Version 6.2

Contents

Introduction.....	3
Compatibility.....	3
Software.....	3
Installation Steps.....	4
1. Importing and loading the taskbar.....	4
2. Running the MX Analysis taskbar installer (Optional).....	7
Multiple Configurations (Groups).....	11
Multi-User mode.....	11
Installation Details.....	12
1. Create recommended folders.....	12
2. Reset file paths.....	12
3. Configure simplified menu.....	12

Introduction

This guide provides instructions on how to install the optional MetaXpress Analysis Taskbar (“MX Taskbar”). The IX Taskbar is an organized collection of tools intended to enhance and/or streamline common tasks and user workflows when reviewing and analyzing images from an ImageXpress high-content screening system.

Compatibility

Software

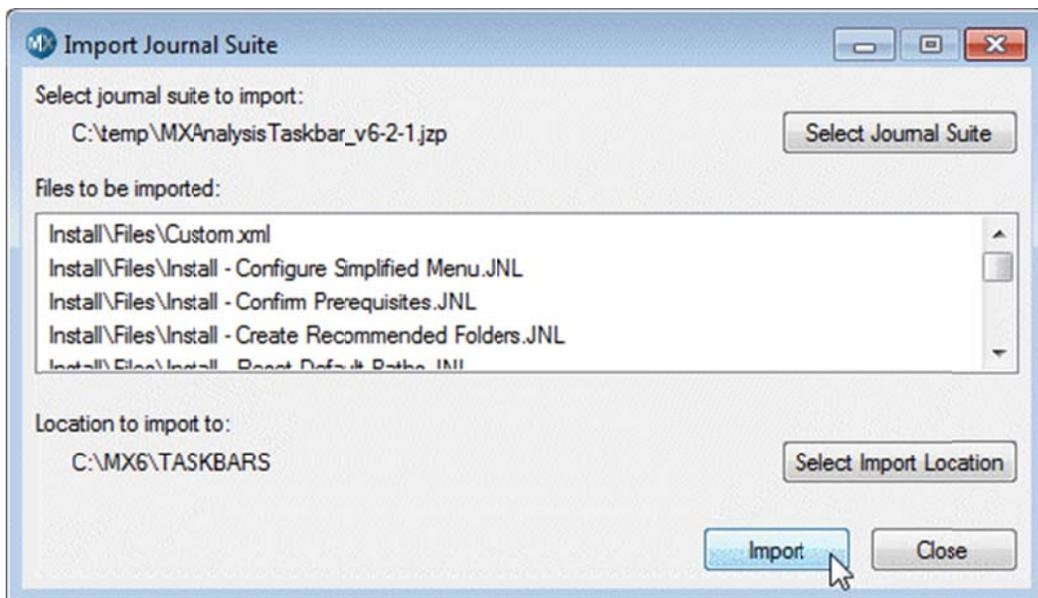
- MetaXpress software version 6.2 and above (offline for analysis)

For ImageXpress acquisition computers, use the IX Taskbar version 6.2, which includes all of the functions in the MX Analysis taskbar.

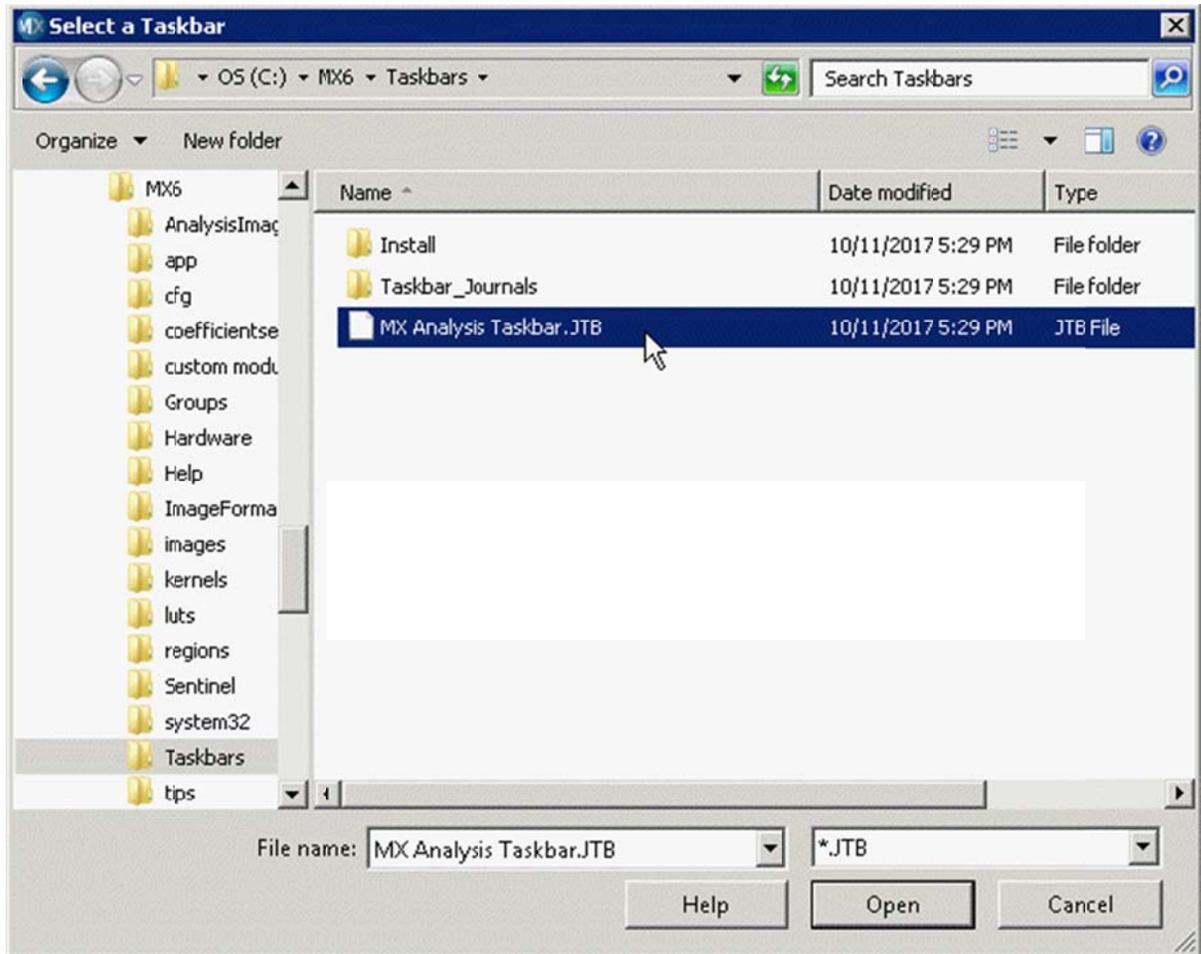
Installation Steps

1. Importing and loading the taskbar

- a. Copy the provided **MXAnalysisTaskbar_v6-2-x.jzp** file to a convenient location on the MetaXpress analysis computer.
- b. Start up the MetaXpress software.
- c. Go to **Journal > Import Journal Suite** (standard menu structure) or **Control > Journal > Import Journal Suite** (simplified menu structure).
- d. Select the **MXAnalysisTaskbar_v6-2-x.jzp** file, and select the “Taskbars” folder within the MetaXpress installation directory (e.g. **C:\MX6\Taskbars**). If the folder does not exist, create it. Click **Import**.



- e. Go to **Journal > Taskbars > Load Taskbar** (standard menu structure) or **Control > Journal > Taskbars > Load Taskbars** (simplified menu structure). Select the **MX Analysis Taskbar.JTB** file in the Taskbars directory (e.g. **C:\MX6\Taskbars**).

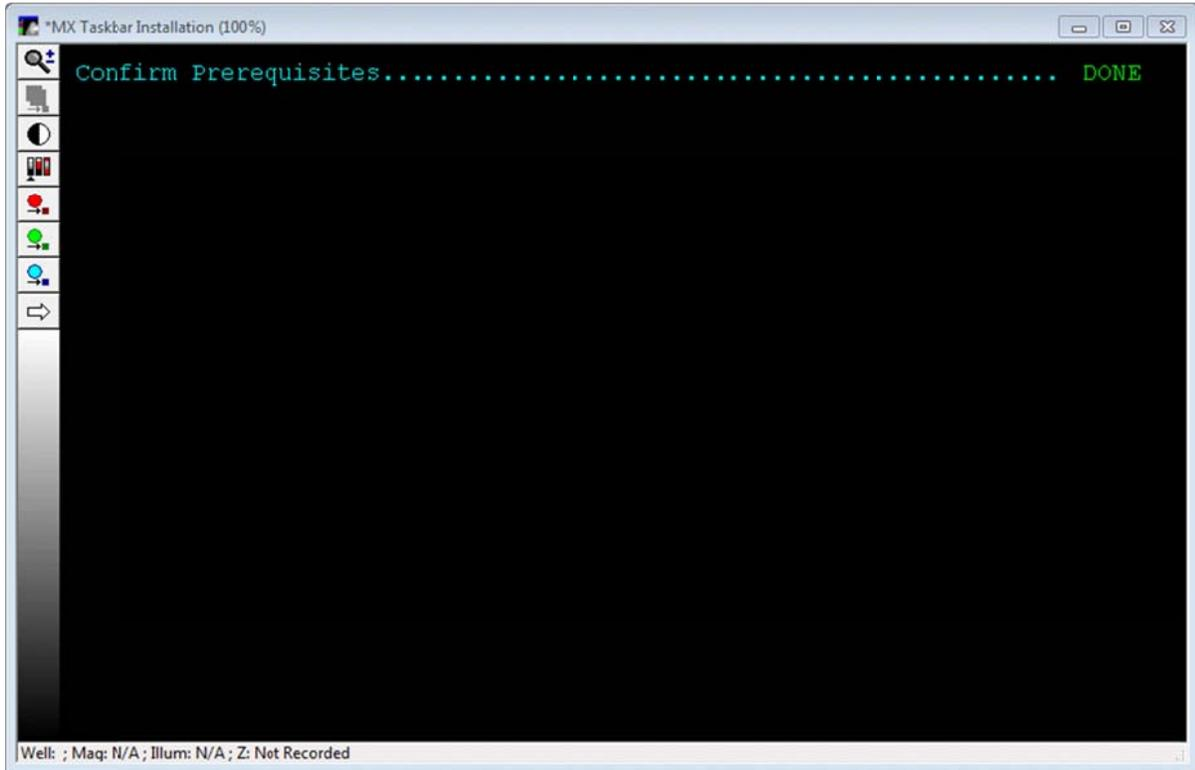


- f. The MX Analysis taskbar appears in the MetaXpress window.

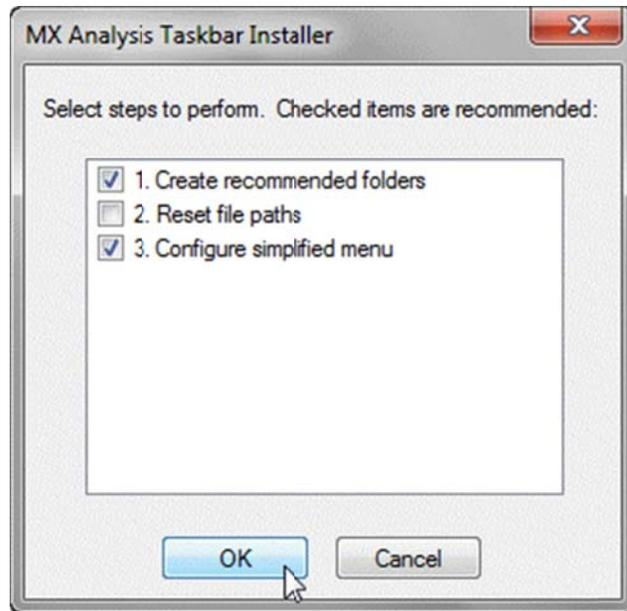


2. Running the MX Analysis taskbar installer (Optional)

- g. On the Main Taskbar, click **Run MX Taskbar Installer (Optional)**.
- h. The MX Taskbar Installation status window appears, confirming if prerequisites have been met:
 - The taskbar files were imported into a valid MetaXpress directory

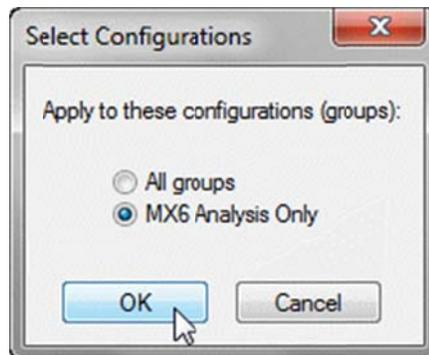


- i. A menu of installation items appears with recommended items already checked. Change the selection as desired, then click **OK**. The individual items are described in detail later in this document.



- j. If either step 2 (Reset file paths) or step 3 (Configure simplified menu) was selected, choose whether to apply the changes to all configurations (also known as groups or profiles) or just the current configuration.

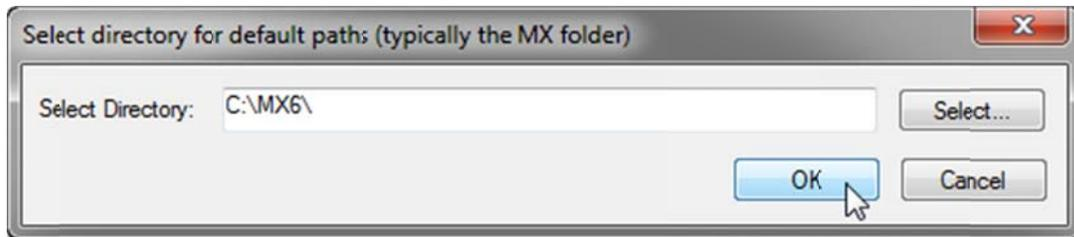
Note: It is uncommon for offline analysis computers to have multiple groups in MetaXpress.



- k. If either step 2 (Reset file paths) or step 3 (Configure simplified menu) was selected, and MetaXpress is running in Multi-User Mode (uncommon), choose whether to apply the changes to all MetaXpress Users (these are different than the database users) or just the current MetaXpress User. If MetaXpress is running in Single-User Mode (default) then this step is skipped.



- l. If step 2 (Reset file paths) was selected, select the base folder to use for the default paths. This is usually either the MetaXpress installation directory (e.g. **C:\MX6**) or a location on the D: drive (e.g. **D:\MX Data**).

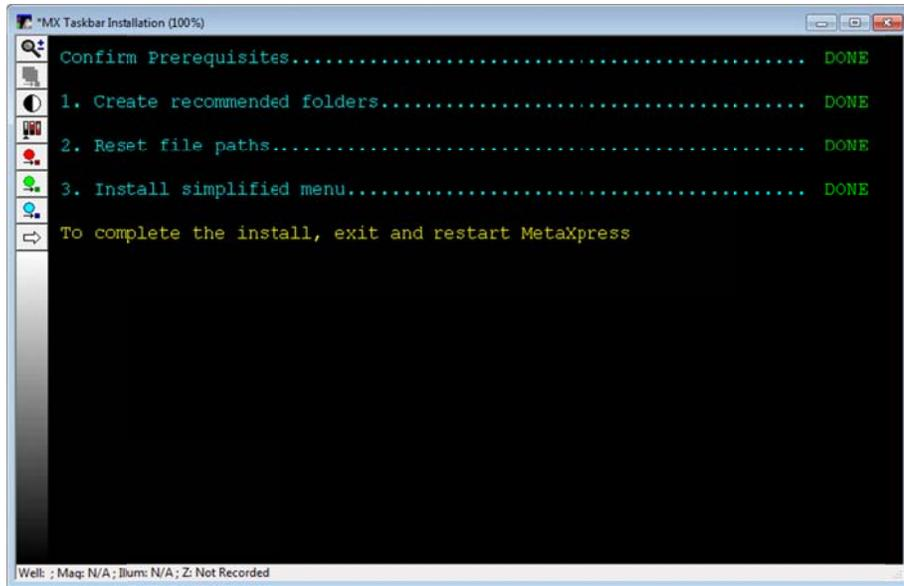


- m. If step 3 (Configure simplified menu) was selected, select whether to install or uninstall the simplified menu.



- n. The installation steps are now performed and the installation status window is updated.

MetaXpress Analysis Taskbar version 6.2 Installation Guide



- o. If step 3 (Configure simplified menu) was selected, exit and restart the MetaXpress software to complete the installation.

Multiple Configurations (Groups)

- p. If there are multiple configurations (multiple MetaXpress program shortcuts), the following step must be repeated for each configuration:
- Load the MX Analysis Taskbar using **Journal > Taskbars > Load Taskbar** (standard menu structure) or **Control > Journal > Taskbars > Load Taskbars** (simplified menu structure)

Multi-User mode

- q. If MetaXpress is run in Multi-User mode (uncommon), the following step must be repeated for each MetaXpress User (this is different than a database user):
- Load the MX Analysis Taskbar using **Journal > Taskbars > Load Taskbar** (standard menu structure) or **Control > Journal > Taskbars > Load Taskbars** (simplified menu structure)

Installation Details

1. Create recommended folders

Step 1 creates the following folders on the computer, if they do not already exist:

C:\assay (required for use of analysis journals)

C:\analysis (required for use of analysis journals)

2. Reset file paths

In the MetaXpress software, the user can use **Edit > Configure Default Paths** to specify the default file paths for image saving, various settings files, regions, etc. However, each path must be specified individually. Step 4 of the taskbar installer allows you to configure all of the default paths in a single step by selecting the base folder to use, which is generally the MetaXpress installation directory (e.g. **C:\MX6**).

NOTE: After installing a new version of the MetaXpress software using the option to copy settings from a previous version, the default paths may be pointing to the previous version installation directory. It is recommended to use the MX Taskbar Installer to update these paths to reduce user confusion.

3. Configure simplified menu

An optional simplified menu structure can be installed to reduce the number of top-level menus. The same features of the software are available in this reorganized menu structure. You can use the Menu Map in the Help menu to help you find the locations of features in the simplified menu structure.

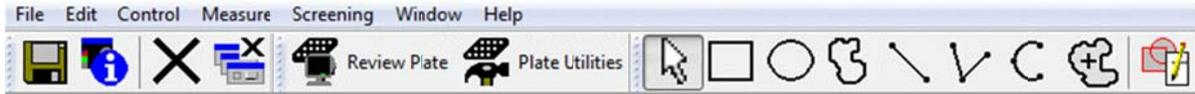
The IX taskbar installer can be used to install or uninstall the simplified menu structure. After either installation or uninstallation, the Metaxpress software must be restarted.

If you already have the simplified menu structure installed, but want to update it to the latest version, choose the install option.

Standard menu structure:



Simplified menu structure:



Right-clicking over the toolbar area will allow you to customize the toolbar selection. The following toolbars are recommended for offline MetaXpress analysis computers:

Basic Tools (enable Show Large Icons)

Screening Tools (enable Show Large Icons and Show Button Text)

Region Tools (enable Show Large Icons, right-click and select Properties to edit Region Tools options)

