



MetaXpress 6 Training

Starting Up and Signing Into MetaXpress



Date Revised 02/03/16 Version B

Index

- [Chapter purpose](#)
- [MetaXpress: Profiles and Shortcuts](#)
- [MetaXpress: Run as Administrator Option](#)
- [MetaXpress: Summary of Startup Steps](#)
 1. [Multi-User Mode: Select User Name](#)
 2. [MetaXpress Splash Screen Appears](#)
 3. [Hardware Initializes](#)
 4. [View License Agreement](#)
 5. [Sign Into the Database](#)
 6. [Size-limited Database Warning](#)
 7. [Select Security Level \(Database Group\)](#)
 8. [StartUp Journal](#)
- [MetaXpress Title Bar](#)
- [Standard or Simplified Menu](#)
- [Toolbars](#)
- [Taskbars](#)
- [MetaXpress Command Line Options](#)
- [Support resources](#)



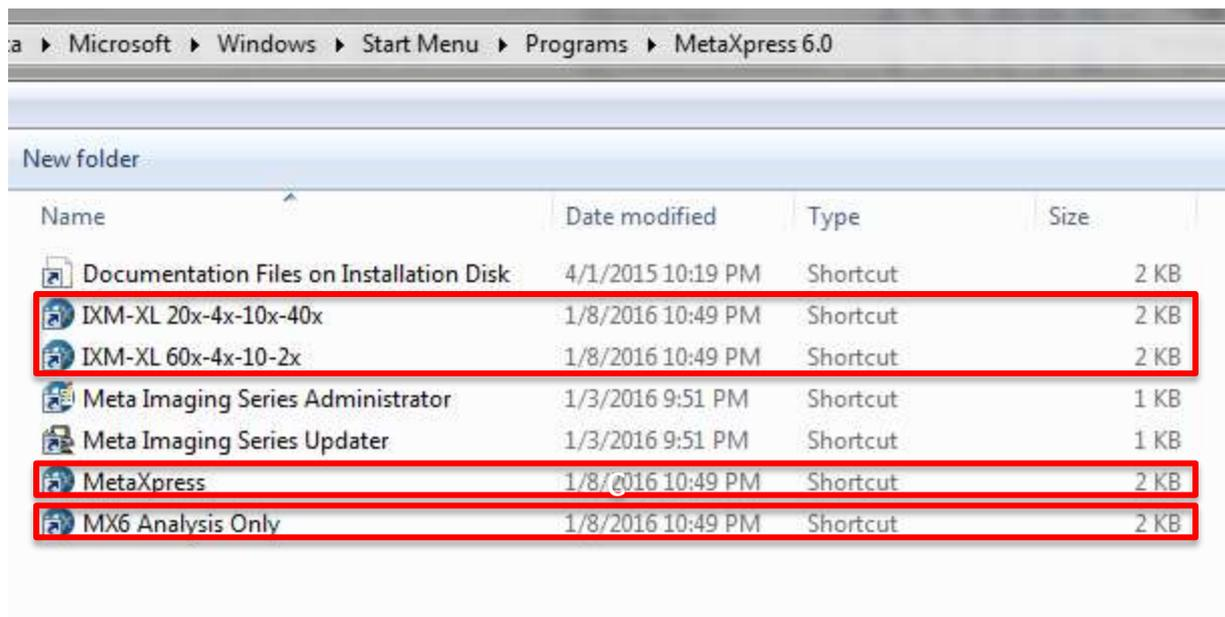
Chapter Purpose

This purpose of this chapter is to explain the different steps and available options when starting up and signing into the MetaXpress software.



MetaXpress: Profiles and Shortcuts

- Some ImageXpress systems have multiple hardware profiles:
 - These are generally named to reflect multiple objective/filter configurations (see example below).
 - Select the shortcut matching your current hardware configuration.
 - You may also see an “offline” or “analysis only” copy – use this to access image analysis functionality without interacting with the IXM hardware.
- Other systems only have a single hardware profile:
 - Use MetaXpress shortcut.



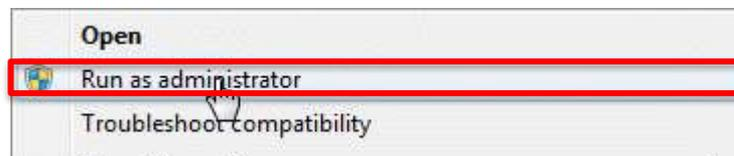
The screenshot shows the Windows Start Menu Programs for MetaXpress 6.0. The path is: Microsoft > Windows > Start Menu > Programs > MetaXpress 6.0. The folder is named "New folder". The contents are as follows:

Name	Date modified	Type	Size
Documentation Files on Installation Disk	4/1/2015 10:19 PM	Shortcut	2 KB
IXM-XL 20x-4x-10x-40x	1/8/2016 10:49 PM	Shortcut	2 KB
IXM-XL 60x-4x-10-2x	1/8/2016 10:49 PM	Shortcut	2 KB
Meta Imaging Series Administrator	1/3/2016 9:51 PM	Shortcut	1 KB
Meta Imaging Series Updater	1/3/2016 9:51 PM	Shortcut	1 KB
MetaXpress	1/8/2016 10:49 PM	Shortcut	2 KB
MX6 Analysis Only	1/8/2016 10:49 PM	Shortcut	2 KB



MetaXpress: Run as Administrator Option

- In some cases it is necessary to run the MetaXpress software “as administrator”:
 - If you want to create a new data source (link to the database).
 - If you are switching between versions of MetaXpress on this computer.
- To run MX as administrator, start the software by right-clicking on the shortcut and selecting the option “Run as administrator.”



- Otherwise, simply double-click on the shortcut.



MetaXpress: Summary of Startup Steps

1. If in Multi-User mode (uncommon): Select user and enter optional password.
2. MetaXpress Splash Screen appears.
3. If connected to an instrument, the hardware initializes.
4. If appropriate, the software license agreement appears.
5. When prompted, sign into the database.
6. If appropriate, the size-limited database warning appears (this will occur if you are running Microsoft SQL Express)..
7. Select the security level for acquisition (hardware profiles only).
8. If a startup journal is selected, it will run.

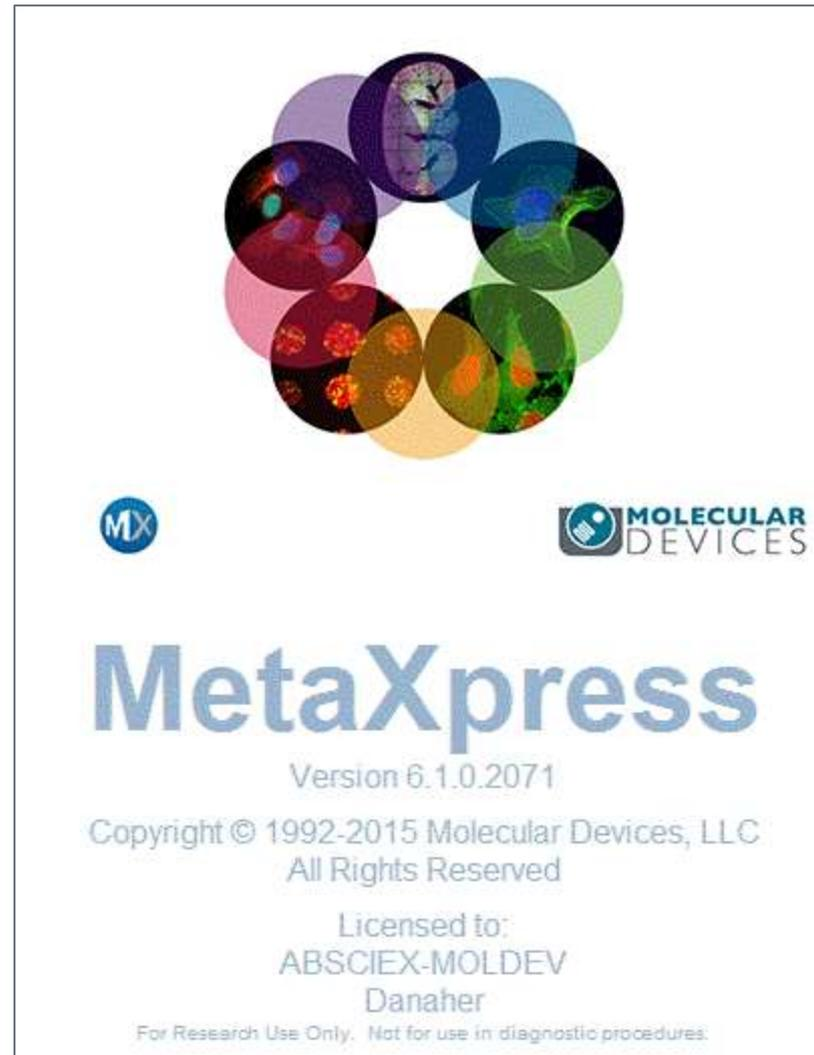


1. Multi-User Mode: Select User Name

- This dialog only appears if MetaXpress is configured in Multi-User mode (uncommon).
- Select your User Name from the dropdown list, then click OK.
- **NOTES:**
 - User Names are not related to database users.
 - Passwords are optional.



2. MetaXpress Splash Screen Appears:



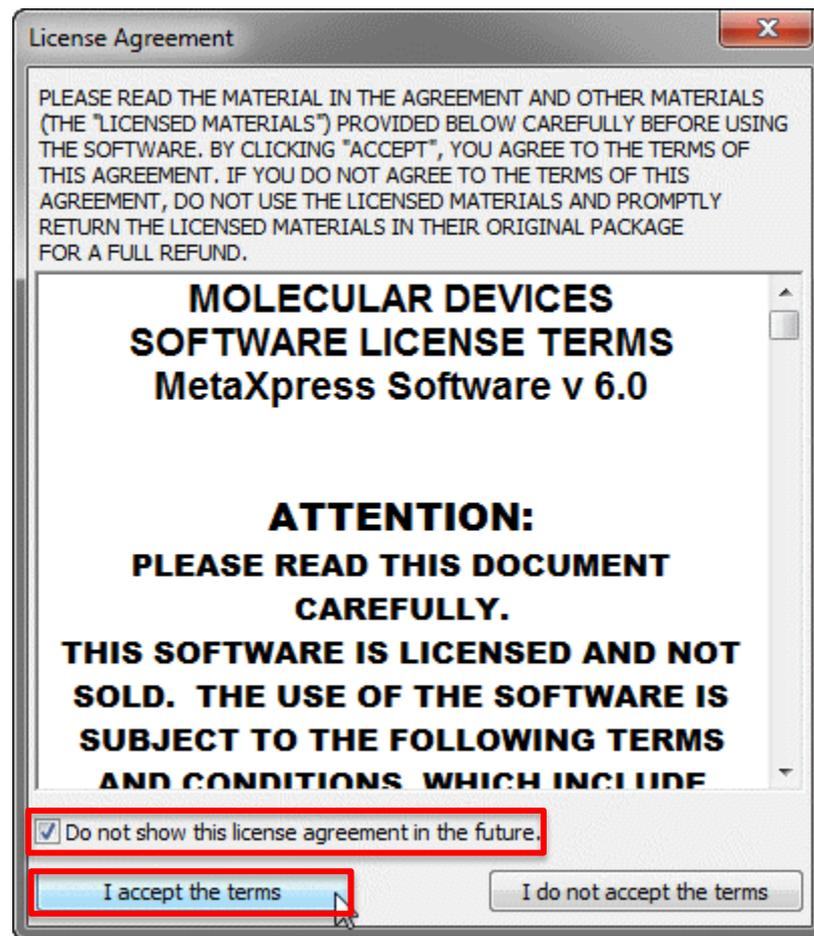
3. Hardware Initializes:

- When the software connects to the IXM instrument for the first time, it homes the various motors.
- Watch for error messages at this point (for example, indicating that the system power is turned off).
- **NOTE: Error messages may become hidden behind the splash screen; to view them, minimize the MetaXpress software window.**



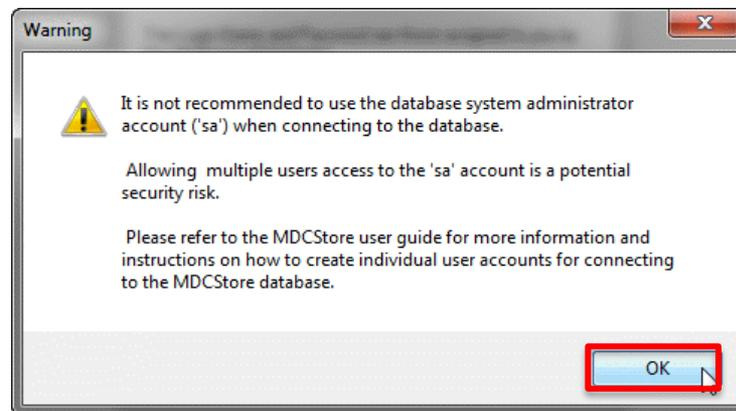
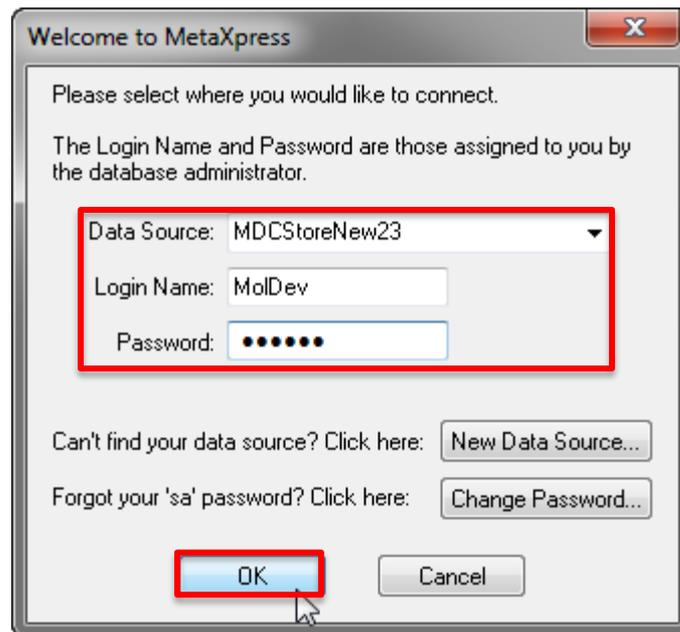
4. View License Agreement

- If this is your first time using the software on this computer, you may see the License Agreement.
- Read, then click on “ I accept the terms” to continue.
- Enable the “Do not show this license agreement in the future” checkbox option to skip this step in the future.



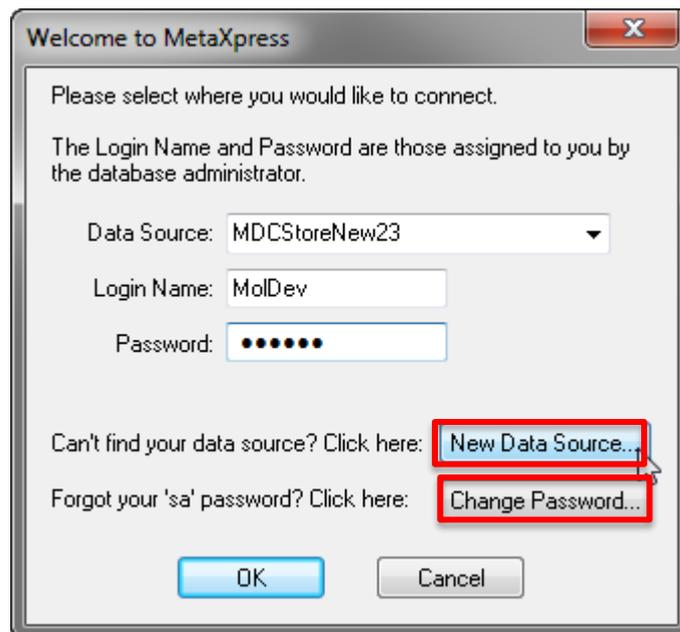
5. Sign Into the Database:

- Select Data Source (link to database) from list.
 - NOTE: Depending on your lab's setup there may be one or many data sources
- Enter database login name and password, then click OK.
- If you use the 'sa' account you will see an additional warning. Click OK.



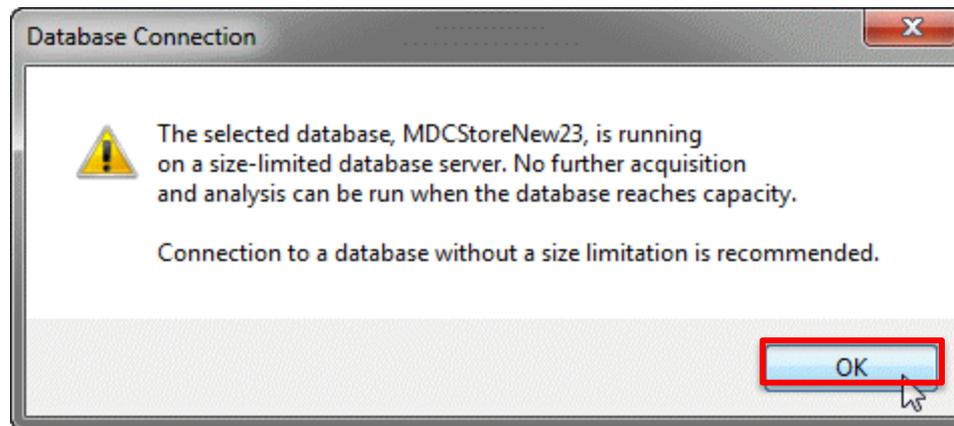
5. Sign Into the Database – Additional Options

- **New Data Source:**
 - Requires launching MetaXpress using **Run as administrator** option.
 - Creates a new link to a database.
- **Change Password**
 - For sa account (MS SQL Server administrative account) ONLY.
 - Requires that the database instance is running on the computer that you are logging into (check with your system administrator to verify).
 - Requires that you are signed into Windows under an administrative account.
 - Does not require old password to be entered.



6. Size-limited Database Warning

- This warning is only displayed if you are using certain database platforms (e.g. SQL Server Express).
- To avoid seeing this warning, upgrade to the full version of the database platform software.
- Click OK to continue.



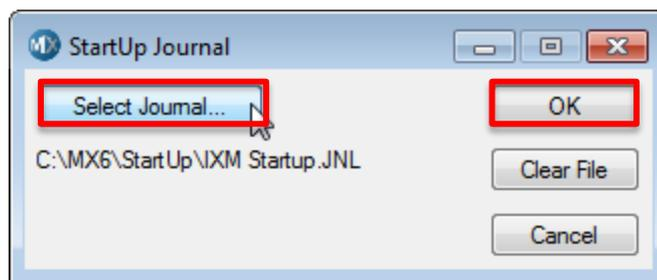
7. Select Security Level (Database Group)

- This dialog is only displayed when launching a hardware configuration.
- Determines which database users will have access to the plates you acquire during this session.
- **NOTE:** Plate security settings can be modified later if necessary.



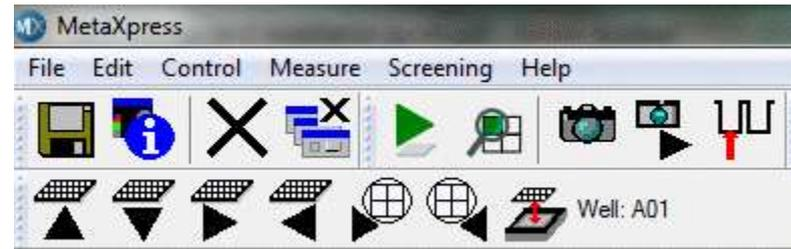
8. StartUp Journal

- Some systems may be configured with an optional StartUp Journal.
- Journal commands start running when the **Select security level** dialog appears.
- Behavior depends on the specific journal.
- A simple example is to load a specific taskbar or specific settings at startup.
- Set StartUp journal in **Journal > Journal Control > Startup Journal** (standard menu) or **Control > Journal > Journal Control > Startup Journal** (simplified menu)

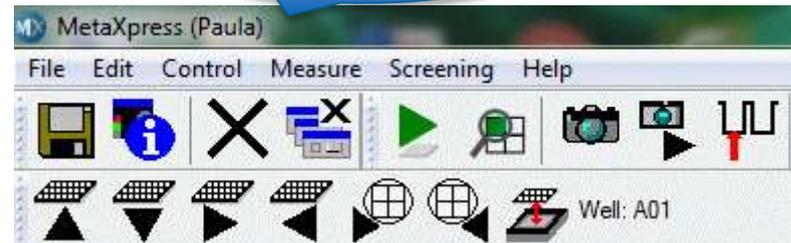


MetaXpress Title Bar

- Single-User Mode
- Multi-User Mode
 - The User Name is shown
- External Control available
 - Used for robot integration
 - External Control will be enabled when the robot software connects to MetaXpress



User Name



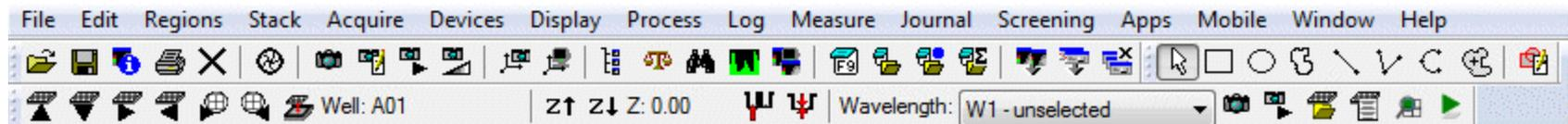
System ID



Standard or Simplified Menu

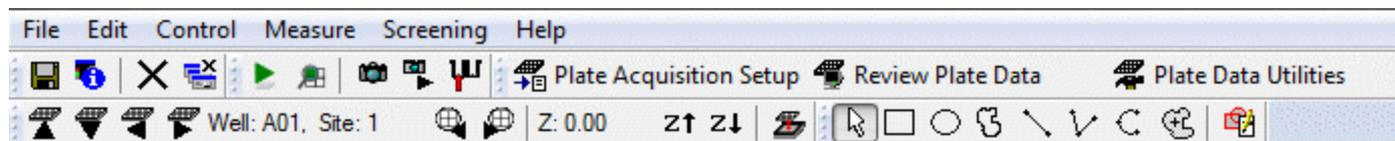
- **Standard Menu:**

- Traditional menu used in earlier versions of the MetaXpress software.



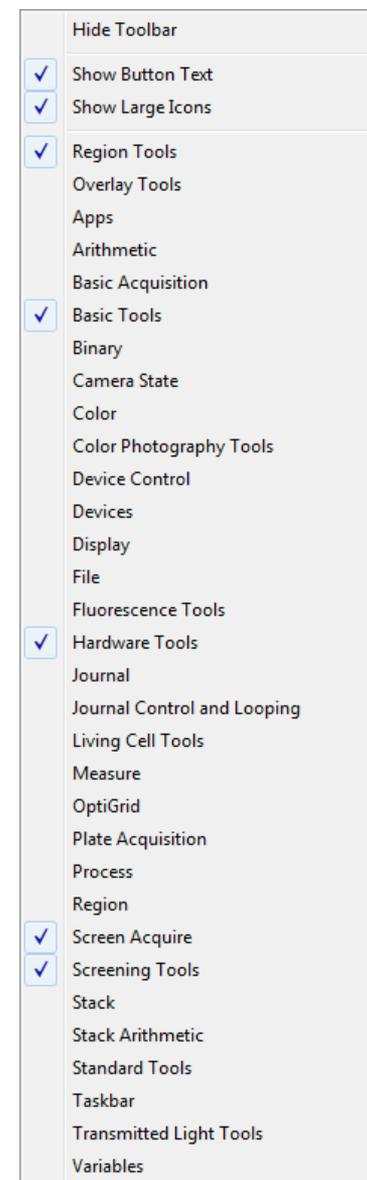
- **Simplified Menu:**

- Introduced with MetaXpress version 6
- Optional.
- Go to **Help > Menu Map** to locate specific commands
- Use the IXM Taskbar Installer to install or uninstall the simplified menu.



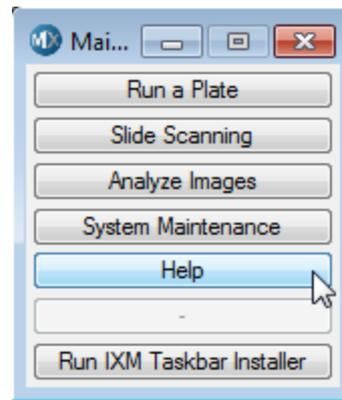
Toolbars

- Toolbars contain links to common functions
- Individual toolbars can be docked horizontally or vertically around any edge of the MetaXpress window.
- Enable or disable specific toolbars by right-clicking on any toolbar or in the toolbar space.
 - Text can be enabled or disabled.
 - Select small or large icons.
- If all toolbars are disabled, go to **Window > Toolbars** or **Control > Window > Toolbars** to access the Toolbar list.
- The Region Tools toolbar can also be enabled from **Regions > Region Tools** or **Measure > Regions > Region Tools**.



Taskbars

- Taskbars are customizable collections of buttons that float on the MetaXpress desktop.
- Taskbar buttons can link to built-in MetaXpress functions, MetaXpress journals, or load other taskbars.
- A standard taskbar is provided with all ImageXpress systems.
- Use the tools in **Journal > Taskbars** or **Control > Journal > Taskbars** to create, edit or load a taskbar.



MetaXpress Command Line Options

- Can be used to reduce the number of steps required for starting the MetaXpress software.
- Available for MetaXpress version 6 and above.
- You can specify any of the following parameters:
 - Data Source Name
 - Login Name
 - Password
 - Database Group
 - MX Profile Name
 - StartUp Journal (runs instead of the StartUp journal selected in the software)
- Use the following format:

```
C:\MX6\mmapp.exe -D<data_source_name> -L<database_login_name>  
-P<database_login_password> -G<database_group_name>  
-J<journal_file_path> <MMAdmin_profile_name>
```

- Enclose the value in quotes if it includes spaces
- Examples:

```
C:\MX6\mmapp.exe -DMDCStore -LMolDev -Pmoldev -GAdministrators  
MetaXpress
```

```
C:\MX6\mmapp.exe -DMDCStore -L"Test User" -P"My Password" -G"Test DB  
Group" "MetaXpress Profile"
```

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



MOLECULAR DEVICES

ADVANCING PROTEIN AND CELL BIOLOGY