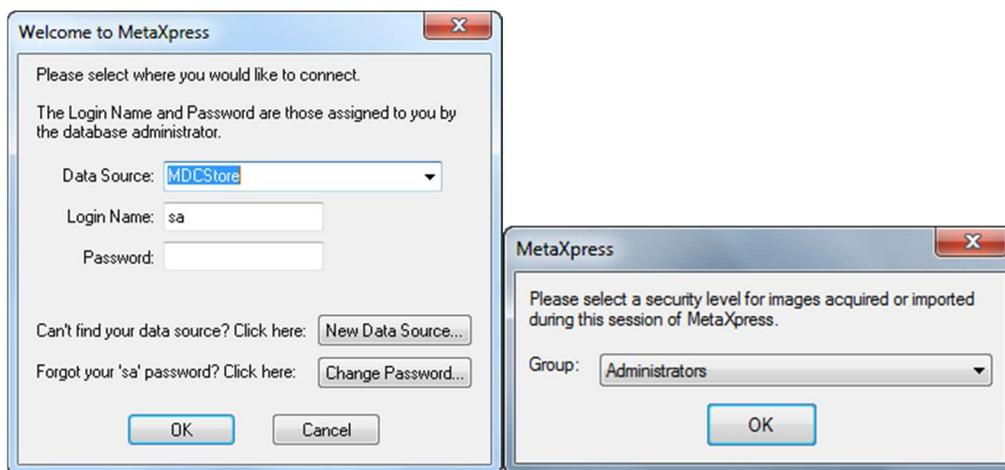


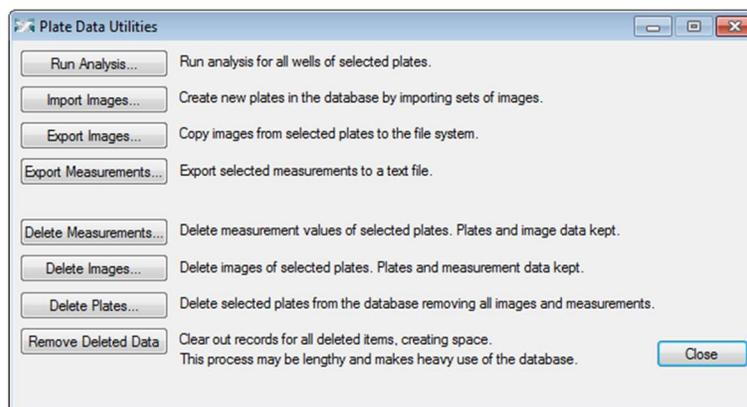
Quick Start Guide – Export Plates (Images) in MetaXpress

Exporting Plates in MetaXpress

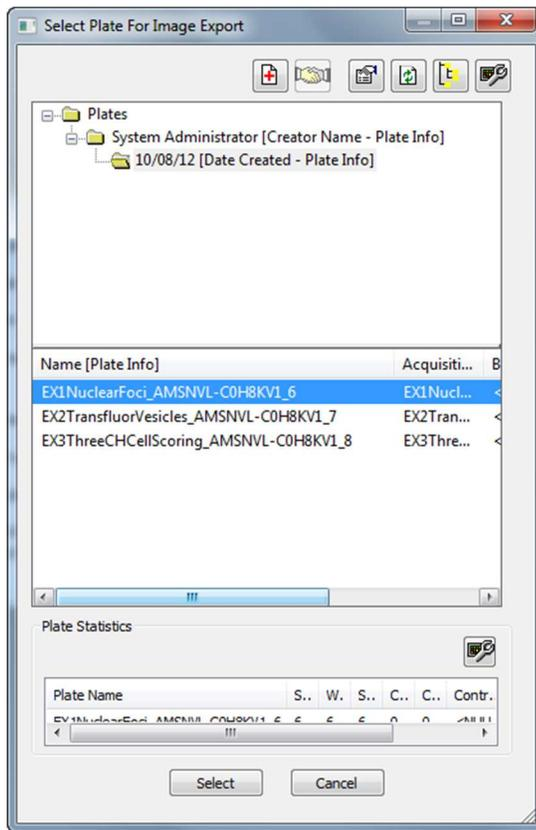
1. Open MetaXpress and log into the database



2. On the main toolbar, click on **Screen → Plate Data Utilities**



3. Click on **Export Images**



4. Highlight the plate(s) you want to export and click **Select**
5. You will then be asked to select a folder in Windows where you would like to save the plate.