



MDCStore™ Xchange Data Conversion Service

Quick Start Guide

5020935 A
August 2012

This document is provided to customers who have purchased Molecular Devices, LLC ("Molecular Devices") equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a license agreement. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Molecular Devices makes no warranties or representations as to the fitness of this equipment for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom.



For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured by Molecular Devices, LLC.
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
Molecular Devices, LLC is ISO 9001 registered.

© 2012 Molecular Devices, LLC.

All rights reserved.

Using the MDCStore Xchange Data Conversion Service

Introduction

The MDCStore™ Xchange Data Conversion Service works together with the installed MDCStore™ High Content Data Management Solution data base to manage image and data conversions.

You can use the built in templates or create custom templates for importing and exporting images and metadata to and from a variety of data sources and image formats.

- Importing supports reading of BMP, JPG, DIB, TIFF, PNG, and GIF formats, as well as OME single and stack image files. You can import image data from HTD, InCell 1000, InCell 2000, Cellomics TIFF, Cellomics DIB, Leica, and BD formats.
- Exporting supports converting image files into BMP (4, 8, 16 bit), JPG (8 bit), DIB (4, 8, 16 bit), Tiff (4, 8, 16 bit), PNG (4, 8, 16 bit), and GIF (8 bit) file formats.



Note: 12 bit images saved as 16 bit Tiffs can be exported only as 16 bit Tiffs.

You can schedule recurring import and export jobs with real-time monitoring of the storage location, and set up email notifications through your SMTP server.

Supported Web Browsers

The MDCStore Xchange Service uses your default web browser to create and schedule import and export jobs. The service supports the following web browsers:

- Microsoft Internet Explorer (version 9 or later)
- Mozilla FireFox
- Apple Safari
- Google Chrome

Installing the Service

The MDCStore Xchange service server and client installers are available on the MetaXpress® Software version 5.0 USB flash drive. For best results, install both the MDCStore Xchange service server and client on your computer. For installation instructions, see the *MetaXpress High Content Image Acquisition and Analysis Software Suite Installation and Update Guide*.

The most recent version of the installer is available on the Molecular Devices support web site. Molecular Devices recommends that you check the support web site and knowledge base for the latest updates to the software and for an updated list of features, supported browsers, and supported operating systems.

To view the Molecular Devices support web site, click the **Support** link at the top of the MDCStore Xchange Service window or go to <http://www.moleculardevices.com/support.html>.

Starting the Service

Before starting the MDCStore Xchange Service, you must set up a data source with the MetaXpress Software or the MDCStore Data Manager. For instructions, see the *MetaXpress High Content Image Acquisition and Analysis Software Suite Installation and Update Guide* or the *MDCStore High Content Data Management Solution Database Schema Installation and Update Guide*.

Running the Background Services

After the MDCStore Xchange Service client and server are installed, the client and server services are started and running in the background. For the MDCStore Xchange Service to function properly, these background services need to be running. If the services have been stopped, you can start them again from the Windows Start menu.

- To start the client service, click **Start > All Programs > Molecular Devices > MDCStore Xchange > Start MDCStore Xchange Client**.
- To start the server service, click **Start > All Programs > Molecular Devices > MDCStore Xchange > Start MDCStore Xchange Server**.

Running the MDCStore Xchange Service

To start the MDCStore Xchange Service in your default web browser, click **Start > All Programs > Molecular Devices > MDCStore Xchange > MDCStore Xchange**.

Logging into your MDCStore Database

In the first MDCStore Xchange service web page, select the **Data Source**, type the **Login Name** and **Password** that you use with the MDCStore Data Manager data base, and then click **Sign In**.

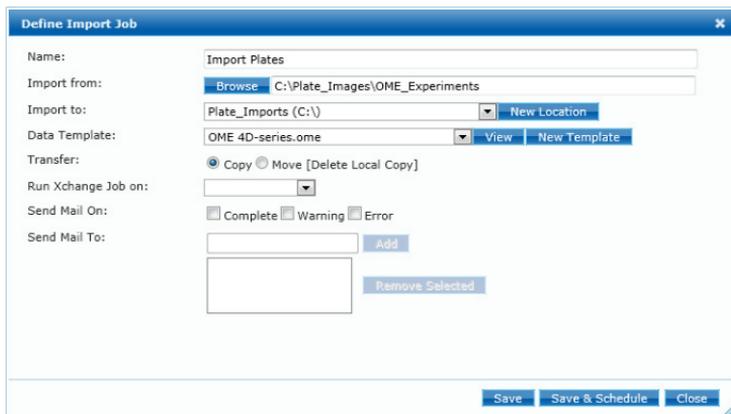
The Login Name and Password are those assigned to you by the database administrator.

Data Source:	<input type="text" value="imageserver1"/>
Login Name:	<input type="text" value="sa"/>
Password:	<input type="password" value="....."/>

[Sign In](#)

Defining an Import Job

1. In the main MDCStore Xchange service window, click **Define Import Job**.
2. In the **Define Import Job** dialog, type a name for the import job.



3. Type the path or browse to the location of the plates to be imported.
4. Define the location to save the imported files.
5. Click the **Data Template** list and select the data template for the import.
The data template defines the layout of the plate data and the conversion method.
6. Click **Copy** to import a copy of the original image. To delete the original image after importing, click **Move**.
7. Click the **Run Xchange Job** on list and select the computer to use for processing the import job.
You can also define email notifications if you have already set up your mail settings. To define your mail settings, click **Settings** in the main MDCStore Xchange service window.
8. Click **Save & Schedule** to open the **Schedule Job** dialog and define when to perform the import. See [Scheduling Import and Export Jobs on page 8](#).

Defining an Export Job

1. In the main MDCStore Xchange service window, click **Define Export Job**.
2. In the **Define Export Job** dialog, type a name for the export job.

The screenshot shows the 'Define Export Job' dialog box with the following details:

- Name:** Export Plates
- Export to:** C:\Plate_Exports (with a 'Browse' button)
- Export from:** (empty dropdown)
- Data Template:** OME Plate Well Site (with 'View' and 'New Template' buttons)
- Data Type:** Images Measurements
- Run Xchange Job on:** IdvWin7Test01 (dropdown)
- Send Mail On:** Complete Warning Error
- Send Mail To:** (empty text field with an 'Add' button)
- Below the 'Send Mail To' field is a list box and a 'Remove Selected' button.
- At the bottom are 'Save', 'Save & Schedule', and 'Close' buttons.

3. Type the path or browse to the location to save the exported files.
4. Click the **Data Template** list and select the data template for the export.
The data template defines the layout of the plate data and the conversion method.
5. Select the **Images** check box to export image data.
6. Select the **Measurements** check box to export measurement data.
7. Click the **Run Xchange Job** on list and select the computer to use for processing the export job.
You can also define email notifications if you have already set up your mail settings. To define your mail settings, click **Settings** in the main MDCStore Xchange service window.
8. Click **Save & Schedule** to open the **Schedule Job** dialog and define when to perform the export. See [Scheduling Import and Export Jobs on page 8](#).

Scheduling Import and Export Jobs

After you define an import or export job, Click **Save & Schedule** to open the **Schedule Job** dialog and define when to perform the job. You can schedule the job to run once or at scheduled intervals.

Scheduling a One-Time Job

1. In the **Schedule Job** dialog, click **One Time**.
2. Click the **Run At** field and select the date and time to run the job.
3. Click **Custom** to define a date range for the source plates. If you want to process all the plates in the source location, then click **All Plates**.
4. Click the **Date From** and **Date To** fields and select the dates and times that define the earliest and latest acquisition dates for the plates to be processed by this job.
5. Select the **Ignore processed plates** check box to prevent processing plates that have already been processed by the MDCStore Xchange service.
6. Click **Save** to schedule the job.

Scheduling a Recurring Job

1. In the **Schedule Job** dialog, click **Recurring**.
2. Click **Add Recurrence**.
3. Click the **Start Time** and **End Time** fields and select the times that define the start and end the job.

Start Time	End Time	Days	
6:00 PM	10:00 PM	<input type="checkbox"/> S <input checked="" type="checkbox"/> M <input checked="" type="checkbox"/> T <input checked="" type="checkbox"/> W <input checked="" type="checkbox"/> Th <input checked="" type="checkbox"/> F <input type="checkbox"/> S	Delete

4. Select the check boxes for the days of the week when you want the job to run.
5. To schedule additional times and days to run this job, repeat [Step 1](#) through [Step 4](#).
6. In **Check for new Plates Every** define the number of minutes between queries for newly acquired plates.
7. Click **Save** to schedule the job.