

GxP Admin Software

Version 3.0.1

User Guide



This document is provided to customers who have purchased 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 non-transferrable license. 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. Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product 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. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

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/patents

Product manufactured by Molecular Devices, LLC. 3860 N. First Street, San Jose, California, 95134, United States of America. Molecular Devices, LLC is ISO 9001 registered. ©2020 Molecular Devices, LLC. All rights reserved.

CE



Contents

Chapter 1: Introduction	5
Primary Data Integrity	5
Administratively Controlled Access	6
Document Workflow	7
Audit Trail	
Password Management	
Chapter 2: Software Installation Introduction	
Backing Up The GxP Admin Database	
Converting Legacy Users	15
Chapter 3: Getting Started	
Logging In To The GxP Admin Portal	
Home - First Steps	20
Chapter 4: Home	21
Change Password	24
Change Date/Time Format	25
Chapter 5: Users	
New Users	
Edit Users	
User Information	
Chapter 6: Roles	
New Roles	
Edit Roles	35
Copy Roles	
Assigned Permissions	
Assigned Projects	
Role Permissions	
Chapter 7: Projects	
New Projects	43
Edit Projects	44
Copy Projects	45
Assign Users to Project Roles	

Project Information	48
Chapter 8: Maintenance	49
Maintenance Settings	50
Maintenance Notifications	51
Maintenance Licenses	52
Maintenance User Import	54
Chapter 9: System Audit Trail	56
System Audit Trail Layout	57
System Audit Trail Filters	58
System Audit Trail Notes	59
System Audit Trail Report	60
Chapter 10: Logging In To The SoftMax Pro Software - GxP Edition	61
Appendix A: Default Roles	63
Appendix B: Audit Events	66
Obtaining Support	71

Chapter 1: Introduction



The SoftMax[®] Pro 7.1.1 GxP Compliance Software Suite controls Molecular Devices[®] microplate readers. This document describes how to use GxP Admin Software whose user interface is the GxP Admin Portal. GxP Admin Software version 3.0.1 stores data documents, audit information, and all user information in a secure Microsoft SQL database. This allows for full control over file administration and permissions of who can do what to files and within the database.

SoftMax Pro Software - GxP edition extends the SoftMax Pro Software - Standard edition into regulated laboratories that work under GMP, GLP, and FDA 21 CFR Part 11 guidelines for secure electronic records. Users must log on to the SoftMax Pro Software - GxP edition with a user name and password. Permissions regulate user actions and user actions trigger events that are recorded to an audit trail for electronic records.



Note: In this user guide, all references to the SoftMax Pro Software - GxP edition use the term SoftMax Pro GxP Software.

This document is specifically for the GxP Admin Software version 3.0.1 and SoftMax Pro GxP Software version 7.1.1.

- The GxP Admin Software version 3.0.1 is not compatible with any SoftMax Pro GxP Software prior to version 7.1.1.
- The SoftMax Pro GxP Software versions prior to version 7.1 must be used with GxP Admin Software version 2.x.
- Ē

Note: The GxP Admin Software has no user interface. The version number for the GxP Admin Software displays in the bottom left corner of the GxP Admin Portal.

See the *SoftMax Pro Data Acquisition and Analysis Software User Guide* and the application help for more information.

Primary Data Integrity

The integrity of raw data are a primary design consideration of SoftMax Pro 7.1.1 GxP Compliance Software Suite. Data that the SoftMax Pro GxP Software acquires are time stamped and traceable to the user who initiates data acquisition. The software saves data in a Microsoft SQL database that cannot be manipulated. Use the System Audit Trail to monitor over 140 events that include, but are not limited to, changes to documents, instrument settings, reduction settings, notes, electronic statements, logins and failed logins, user activities, and file creation, modification, and deletion.

Documents that the SoftMax Pro GxP Software creates are strictly bound to features that support FDA 21 CFR Part 11. As a result, these documents cannot be opened or modified in the SoftMax Pro Software - Standard edition to ensure the integrity of the FDA 21 CFR Part 11 compliant data.

Ē

Note: No software by itself can be FDA 21 CFR Part 11 Compliant. Use the SoftMax Pro 7.1.1 GxP Compliance Software Suite features to demonstrate compliance with these regulations.

Administratively Controlled Access

The GxP Admin Portal Software is the interactive user interface for the GxP Admin Software. Use the GxP Admin Portal to create users. You grant each user Access permissions to allow them to use the software applications to which they need access. You then assign each user with the SoftMax Pro Access permission to Projects that limit their access to documents. For each Project, you assign the user to a Role that allows them to perform the tasks for which they are responsible.

The computer that runs the SoftMax Pro GxP Software must be able to access the server that contains the GxP Admin Software to enforce the features that are under administrative control.

Application Access Permissions

Use the three Access permissions to restrict each user's access to the software applications.

- **GxP Admin** Users with this Access permission have access to all GxP Admin Portal functionality. GxP Admin Access permission allows management of the users of both software applications including workflow permissions and password resets. GxP Admin users manage application setting such as password strength requirements and event notifications. Users with this Access permission have the System Audit Trail Access permission by default and can view the System Audit Trail in the GxP Admin Portal. These users cannot log on to the SoftMax Pro GxP Software unless you also grant them the SoftMax Pro Access permission.
- System Audit Trail Users with this Access permission can view the System Audit Trail in the GxP Admin Portal. These users can confirm that the GxP Admin Software meets compliance requirements but cannot change user settings, cannot change application settings, and cannot log on to the SoftMax Pro GxP Software unless you also grant them the GxP Admin Access permission and/or the SoftMax Pro Access permission.
- **SoftMax Pro** Users with this Access permission also require a user license to log on to the SoftMax Pro GxP Software. These users are granted permission to access the SoftMax Pro GxP Software based upon their assignment to Projects and their Role in each Project. These users cannot log in to the GxP Admin Portal unless you also grant them the System Audit Trail Access permission and/or the GxP Admin Access permission.

Projects - Document Access Restrictions

Each user with the SoftMax Pro Access permission must be a member of at least one Project. Use Projects to limit which documents the user can access. Each user can be a member of multiple Projects.

Roles - Workflow Permissions

Each user with the SoftMax Pro Access permission can have one Role in each Project. Each user can be a member of multiple Projects and can have a different Role in each Project. Use Roles to restrict the user's responsibilities in the SoftMax Pro GxP Software.

Document Workflow

The document life-cycle in a compliant lab consists of multiple different development and review steps and requires a system to indicate the document state. The status system allows a Project team to track and approve documents while the document moves through the pipeline of development, review, release and usage in a controlled environment.

There are two types of documents:

- Protocols contain instrument settings, notes, statements, group settings, graph settings, and so on. Protocols do not contain data.
- Data documents contain everything from the protocol plus the data that the instrument collects and the results of the formula calculations you set up in the protocol.

Document Statuses

The SoftMax Pro 7.1.1 GxP Compliance Software Suite installation includes three default Roles; Scientist, Lab Technician, and Lab Manager. The default permissions for these Roles correspond to the document statuses that are present in the SoftMax Pro GxP Software. You can modify the default Role permissions, if needed.

- In Work
- Review Pending
- Reviewed
- Released
- In Progress
- Approval Pending
- Approved
- Outdated
- Canceled

The following document workflow corresponds to the default Roles. Your laboratory can define Roles and document workflows that meet your needs.



In Work Document Status - Scientist Role

The scientist opens and saves a new data document. They define the instrument settings in Plate sections and define data analysis settings. The software assigns a new document the status In Work. The scientist adds statements to the document to provide communication to the other users responsible for the document. The scientist signs the first statement to prevent further edits and sets the document status to Review Pending.

Review Pending Document Status - Lab Manager Role

The lab manager reviews the data document with the status Review Pending, signs the second statement, and sets the document status to Reviewed.

Reviewed Document Status - Release Protocol - Lab Manager Role

If the data document with the status Reviewed is correct, the lab manager then releases the document as a protocol. The software saves the protocol with the status Released for reuse. If the document has a newer iteration or is incorrect, the lab manager can set a document that has the status Released to the Outdated status.

In Progress Document Status - Lab Technician Role

The lab technician opens a protocol with the Released status and the software prompts the lab technician to save the protocol as a data document before they can run the experiments. Because the Lab Technician Role has the Generate Compliance Data permission, the software sets the data document status to In Progress and removes the two signed statements. This allows the lab technician to run the experiments and generate the compliance data.

When the lab technician finishes their work, they sign a statement to prevent further changes and set the document status to Approval Pending.

Approval Pending Document Status - Lab Manager Role

The lab manager confirms that the document with the Approval Pending status meets regulatory requirements, signs a statement, and sets the document status to Approved.

Canceled Document Status

Users with appropriate permissions can set the document status to Canceled at the applicable points in the document workflow.

Audit Trail

The GxP Admin Software and the SoftMax Pro GxP Software generate audit trails that record the events that user actions trigger. For each event instance, the audit trail records the date and time, the username of the person who was logged on, and the event action. Users cannot alter the audit trail, nor can they edit or delete audit trail entries.

GxP Admin Software Events

The GxP Admin Software saves the audit trail to the database and you view the System Audit Trail in the GxP Admin Portal.

For a list of the events that trigger a System Audit Trail entry, see Audit Events on page 66.

SoftMax Pro GxP Software Events

In addition to the GxP Admin Software System Audit Trail, the SoftMax Pro GxP Software records document-specific events to an audit trail for the document. You view the document-specific audit trail in the SoftMax Pro GxP Software.

- For protocols, the SoftMax Pro GxP Software does not retain an audit trail. The GxP Admin Software System Audit Trail records applicable protocol event information.
- For data documents, the SoftMax Pro GxP Software audit trail records events that pertain to the document that is active in the Workspace and saves the document-specific events with the data document. Additionally, the document-specific audit trail includes general activities such as log in and log out when a document is open during those events.

The audit trail for each Plate section in a data document includes the following information:

- The name of the user who logged in when the read began.
- The instrument type used to do the read and the instrument ROM version number. For the SpectraMax[®] i3x Multi-Mode Microplate Reader and the SpectraMax[®] Paradigm[®] Multi-Mode Microplate Reader, the audit trail also includes the serial number of the detection cartridge.

Password Management

During installation, you create the first user that you will use to login to the software. This user has the GxP Admin Access permission and the System Audit Trail Access permission. The multi computer setup installation supports the creation of both Custom Credentials users and the integration with Windows Credentials, Active Directory (LDAP) users.

- **Custom Credentials**: For users you create with the Custom Credentials login mode, the software creates a randomly generated alphanumeric string of characters to be their password. This password grants users with the appropriate Access permissions access to the GxP Admin Portal Software and the SoftMax Pro GxP Software. Both software applications require Custom Credentials login mode users to change their password upon initial login. Both software applications enable Custom Credentials login mode users to change their password as often as they want. The GxP Admin Portal has settings to manage password strength and password aging to require Custom Credentials login mode users to change their password at the interval you specify.
- Windows Credentials: For users you create with the Windows Credentials login mode, password management is maintained through the Active Directory. Windows Credentials login mode users cannot change their password in the GxP Admin Portal Software or the SoftMax Pro GxP Software. The password maintenance settings in the GxP Admin Portal do not affect Windows Credentials login mode users. Windows Credential login mode users with the SoftMax Pro Access permission must have a license and must be a member of at least one Project. Their activity within the SoftMax Pro GxP Software is restricted by their Role permissions. Windows Credentials login mode is available when you use the multi computer setup installation.

Chapter 2: Software Installation Introduction



The SoftMax Pro 7.1.1 GxP Compliance Software Suite scales with the size of your laboratory environment. Depending on the size of your implementation, you can install the interdependent SoftMax Pro 7.1.1 GxP Compliance Software Suite components on multiple networked computers. For the smallest of implementations, it is possible to install the components on a single computer. You should consult with your Molecular Devices representative and your network administrators to plan and scale your implementation according to your requirements.

SoftMax Pro 7.1.1 GxP Compliance Software Suite components:

- GxP Admin Software: The GxP Admin Software (server) is the database and security back-end application that you install on a dedicated, centrally located server with a static IP address. This server must be accessible to all computers that run the GxP Admin Portal Software and the SoftMax Pro GxP Software. The GxP Admin Software does not have a user interface.
- **GxP Admin Portal Software**: The GxP Admin Portal is the user interface that you use to interact with the GxP Admin Software. Install the GxP Admin Portal Software on a computer or computers that have access to the dedicated server on which you install the GxP Admin Software.
- **SoftMax Pro GxP Software**: The SoftMax Pro GxP Software (client) controls the microplate reader for data acquisition and statistical data analysis. These computers must have access to the dedicated server on which you install the GxP Admin Software.

Note: In this document, all references to the SoftMax Pro Software - GxP edition use the term SoftMax Pro GxP Software.

See the following documents for installation instructions:

- SoftMax Pro Data Acquisition and Analysis Software GxP Edition Installation Guide for the Multi Computer Setup
- SoftMax Pro Data Acquisition and Analysis Software GxP Edition Installation Guide for the Single Computer Setup

Backing Up The GxP Admin Database

For the single computer setup, the SoftMax Pro 7.1.1 GxP Single Computer Setup Installation Wizard allows you to install the GxP Admin Backup Tool. If you do not have standard Microsoft SQL backup tools, you can use this tool to schedule recurring database backups and to do a backup on an as needed basis.

Note: The GxP Admin Backup Tool is not applicable for the multi computer setup because it is assumed that your network has standard backup procedures in place.

If a backup occurs while the system is in use, there is the possibility that the backup audit trail information could be inconsistent. For example, if a user is actively creating new plates when the backup starts, the backup may save the audit trail entry before the changes to the document are saved. The user continues to work, and the active database is fully consistent. However, the backup contains the audit trail information for the new plates and does not contain the document version with the latest plate added. This is only relevant if you restore the database from such a backup.



Note: Backing up to your Windows user folders, such

C:\Users\<CurrentUser>\Documents is not supported and will result in a backup error. Please choose C:\Users\Public\Documents, or other location not in the Windows Users subdirectory.

Backing Up The Database

The GxP Admin Backup Tool creates a report that documents the relevant information in the backup destination folder.

To back up the GxP Admin Software database:

- 1. Click Or locate and double-click **GxPAdminBackupTool.exe** to display the GxP Admin Backup Tool dialog.
- 2. In the **Backup** area, in the **Destination** field either enter the path to the location to where

you want to back up the database or click 🛄 to display the Browse For Folder dialog.

- 3. In the Browse For Folder dialog, select the destination location and click **OK**.
- 4. Click Backup Now.

Note: If you logged into the Windows operating system as a user that has administrator privileges to the computer, the GxP Admin Backup Tool displays "Could not backup database" if you select to backup to a folder to which you do not have write permissions such as the Documents folder.

- 5. The **History** area displays the operation, status, operator, file name (file name is the date and time the backup is created and includes the path to the file), start time, duration, and size of the backup file. The GxP Admin Backup Tool creates a ReportDocx.pdf file in the backup folder for compliance purposes.
- 6. To manage the backups that display in the History area, click **Settings** to display the Keep Backups dialog.
- 7. In the Keep the Last ___ Backups field, enter the number of backups to save and click OK.
 - Note: If you set this to 0 (zero), no backups will be saved and all existing backups will be deleted.

Creating a Backup Schedule

Use the GxP Admin Backup Tool and Windows Scheduler to schedule regular database backups.

To create a database backup schedule:

- 1. Click Or locate and double-click **GxPAdminBackupTool.exe** to display the GxP Admin Backup Tool dialog.
- 2. In the **Backup** area, click **Create Schedule** to display the Create Schedule Information dialog.
- 3. Click **Open Windows Scheduler** to display the Task Scheduler dialog.
- 4. On the left, select **Task Scheduler Library**.
- 5. In the Task Scheduler Library, select the GxPDatabaseBackupTask.
- 6. On the right in the Actions area, click Properties to display the Properties dialog.
- 7. Select the **Triggers** tab.

- 8. Use the **New**, **Edit**, and **Delete** buttons to define when to schedule the backup task to trigger.
- 9. In the Properties dialog, select the General tab.
- 10. In the **Security Options** area, note the name that appears under When Running The Task, Use The Following User Account. Click **Change User or Group** to select a different user account, if needed.
- 11. Choose an option:
 - Select **Run Only When User Is Logged On** to have the backup schedule task run when a user is logged on.
 - Select **Run Whether User Is Logged On Or Not** (recommended) to have the backup schedule task run without the need for the user to log on. The user you select in the previous step needs to have the LocalSecurityPolicy assignment "Log on as batch job" Windows permission.
- 12. In the Properties dialog, click **OK**. If you select Run Whether User Is Logged On or Not, enter the user's password.
- 13. Close the Task Scheduler dialog.

Restoring From Backup

Note: When you restore the database from a backup, the backup files replace all of the files in the GxP Admin Software database.

To restore the GxP Admin Software database from a backup:

- 1. Click Solution or locate and double-click **GxPAdminBackupTool.exe** to display the GxP Admin Backup Tool dialog.
- 2. In the **Restore** area, in the **Restore From** field, either enter the path to the location of the backup file or click to display the Browse For Folder dialog.
- 3. In the Browse For Folder dialog, select the backup file and click **OK**.
- 4. Click Restore Now.
- 5. Confirm that no other user is logged into the GxP Admin Portal Software or the SoftMax Pro GxP Software.

Note: Whenever you restore the database, you must confirm that no users are using the SoftMax Pro GxP Software or the GxP Admin Portal Software. Restoring while users are logged in may lead to inconsistent data.

- 6. Click **OK** in the message dialog. The History area displays the operation, status, operator, file name, start time, duration, and size of the backup file.
- 7. After a restore operation, check the recent audit entries for inconsistencies and make sure to document any inconsistencies accordingly.

Converting Legacy Users

The GxP Admin Software version 3.x is significantly different from the previous versions. The database structure is different and additional security permissions have been implemented. Therefore, if you want to use the users from GxP Admin Software version 2.x, you must convert legacy user files to an updated format and then import the legacy users into the GxP Admin Software version 3.x database.

Note: If use the multi computer setup and you intend have users use your company's Active Directory database you will use the GxP Admin Portal to create new Windows Credentials login mode users. GxP Admin Software version 2.x did not support Windows Active Directory and users you convert are created as Custom Credential users in the GxP Admin Software 3.x software. The single computer setup does not support the Windows Credentials login mode.

The GxP Admin EDB Converter Tool is a separate software application that converts the version 2.x user files that have an .edb file extension to the version 3.x user files that have an .edbexp file extension.

Note: After you convert the user files, you use the GxP Admin Portal to import usernames, passwords, and the GxP Admin Access permission into the GxP Admin Software version 3.x database. You must then assign each user with the SoftMax Pro Access permission a license from the User Administration page and to a Project from the Project Administration page.

To prepare to convert GxP Admin Software version 2.x user files that have the .edb file extension:

- You can run the GxP Admin EDB Converter Tool on the computer on which you install the SoftMax Pro 7.1.1 GxP Compliance Software Suite if the computer has a network connection to the computers that contain the .edb files.
- You can use a flash drive to copy the .edb files to any location on the computer on which you install the SoftMax Pro 7.1.1 GxP Compliance Software Suite.
- You can use a flash drive to move the GxP EDB Converter folder to the computer on which the .edb files are located.
- You can install the GxP Admin EDB Converter Tool on other computers.

Install the GxP Admin EDB Converter Tool

To install the GxP Admin EDB Converter Tool:

- 1. Insert the GxP Admin Software DVD in the computer DVD drive or use alternative media. Locate and double-click EDBConverter<n.n>Setup.exe file to start the install.
- 2. On the Welcome page, click Next.
- 3. On the License Agreement page:
 - a. Select the I Accept the Agreement check box.
 - b. Click Next.
- 4. On the Destination Folder page and the pages that follow, click **Next** to accept the default settings or you can read the instructions and do a custom installation until the Installation Complete page displays.
- 5. On the Installation Complete page, click **Finish**.

The GxP Admin EDB Converter Tool icon Appears on the computer desktop.

Converting Legacy User Files

The GxP Admin EDB Converter Tool is a separate software application that you run on a computer that has access to your legacy GxP Admin Software user files. The GxP Admin Software version 2.x user files have an .edb file extension and the GxP Admin Software version 3.x user files have an .edbexp file extension.

8 Converter	_		\times
Convert Options			
C Single File All Files in a Folder			
File or Folder Name		_	
Include all Subfolders			
Processing			
Log		_	
Сору			
	Convert	Can	el

Converting a Single File

To convert one .edb file to an .edbexp file:

- 1. Click or locate and double-click **ConverterTool.exe** to display the Converter dialog.
- 2. Select the Single File Convert Option.
- 3. To the right of the File or Folder Name field, click to display the Open dialog.
- 4. Navigate to the .edb file and click **Open** to display the path to the file in the File or Folder Name field. Note this location because the software places the new .edbexp file here.
- 5. Click **Convert**. The Log field displays the progress of the conversion.
- 6. The software places the .edbexp file in the same folder as the .edb file. Copy the .edbexp file to a flash drive or place it on a network drive that is accessible to the computer running the GxP Admin Portal Software. See Maintenance User Import on page 54.

Converting Multiple Files

To convert multiple .edb files to .edbexp files:

- 1. Click or locate and double-click **ConverterTool.exe** to display the Converter dialog.
- 2. Select the All Files In a Folder Convert Option.
- 3. To the right of the File or Folder name field, click to display the Browse For Folder dialog.
- 4. Navigate to the folder that contains the .edb files and click **OK** to display the path to the folder in the File or Folder Name field.
- 5. Select the **Include All Subfolders** check box to convert all .edb files in the subfolders of the folder you select.
- 6. Click **Convert**. The Log field displays the progress of the conversion. The EDB Converter Tool creates an .edbexp file for each .edb file in the same location as the .edb file.
- 7. Copy the .edbexp files to a flash drive or place them on a network drive that is accessible to the computer running the GxP Admin Portal Software. See Maintenance User Import on page 54.

Chapter 3: Getting Started



Now that you have installed the GxP Admin Portal, click On the computer desktop to display the GxP Admin Portal Login page. After you login, you add the licenses to assign to the users to whom you grant the SoftMax Pro Access permission and then add users.

Logging In To The GxP Admin Portal

Use the Login page to access the GxP Admin Software. Users with the GxP Admin Access permission and/or the System Audit Trail Access permission can use the GxP Admin Portal to login to the GxP Admin Software.

C GxP Server Setup	GxP Admin Portal	
GxP Server Setup	GxP Admin Portal	
	Login	$\overline{)}$

Setting Up The GxP Server

The first time you use the GxP Admin Software you need to connect the GxP Admin Portal to the GxP Admin Software database. You do this one time for each GxP Admin Portal you use to administer the GxP Admin Software.

To connect the GxP Admin Portal to the GxP Admin Software:

- 1. On the top left of the Login page, click **GxP Server Setup** to display the GxP Server Hostname page.
- 2. In the Server Hostname field:
 - For the multi computer setup installation, enter either the hostname or IP address of the computer on which you install the GxP Admin Software. The GxP Admin Software computer is NOT the external Microsoft SQL database server if you configure the GxP Admin Software to use an external Microsoft SQL database.
 - For the single computer setup installation, enter **localhost**.
- 3. Click **Continue** to return to the Login page.

Logging In For The First Time

During software installation, you create the first user that you use to login to the software. This user has the GxP Admin Access permission and the System Audit Trail Access permission. Upon first login, this user must change their password.

Note: You get five attempts to log in for the first time. If you forgot the username and password you entered during installation, you need to reinstall the software.

The software assigns all users you create with the Custom Credentials login mode a randomly generated alphanumeric string as their initial password. The software prompts users with the Custom Credentials login mode to change their password upon initial login. Default password strength settings require you to create a password that contains at least ten characters, both uppercase and lowercase letters, at least one digit, and at least one special character. You can change the default password strength settings on the Maintenance page. See Maintenance Settings on page 50.

Users you create with the Windows Credentials login mode use the password associated with their Active Directory user account. Windows Credentials login mode users cannot change their password in the GxP Admin Portal and the software does not prompt these users to change their passwords upon initial login. Windows Credentials login mode user passwords are managed via your company's Active Directory and the password management settings in the GxP Admin Portal are ignored.

To login for the first time:

- 1. In the User Name field enter your user name.
- 2. In the **Password** field, enter your password.
- 3. Click Login.

For Windows Credentials login mode users, the Home page displays. For Custom Credentials login mode users, the Change Password dialog displays.

- 4. For Custom Credentials login mode users, in the **New Password** field, enter your new password with at least ten characters, both uppercase and lowercase letters, at least one digit, and at least one special character.
- 5. In the **Repeat New Password** field, reenter the password.
- 6. Click **Change** to display the Home page. See Home on page 21.

Home - First Steps

When you login to the GxP Admin Portal for the first time, the Home page displays a list of tasks you should do as the first steps.



The general sequence of steps to get started follows the sequence of the items in the First Steps list.

- Add Licenses: Each user with the SoftMax Pro Access permission requires a license to login to the SoftMax Pro GxP Software. See Maintenance Licenses on page 52.
- Import Legacy Users: If you use GxP Admin Software version 2.x, this option allows you to import your existing users into GxP Admin Software version 3.x. See Maintenance User Import on page 54.
- **Create Users**: Use to create the users for both the GxP Admin Software and the SoftMax Pro GxP Software. See Users on page 26.
- **Change Settings**: Use to define settings for password strength and administrator contact information. See Maintenance Settings on page 50.
- **Manage Roles**: Use Roles to restrict what actions users with the SoftMax Pro Access permission can do in the SoftMax Pro GxP Software. See Roles on page 33.
- **Create Projects**: Use Projects to restrict which documents users with the SoftMax Pro Access permission can access in the SoftMax Pro GxP Software. See Projects on page 42.

To permanently remove the First Steps check list, click Get Started.

Note: After you click Get Started you cannot display the First Steps check list again.

Chapter 4: Home



The Home page displays when you login and provides shortcuts to most workflows along with other useful information. The top right of most pages displays the Help icon and the name of the user that logged in. The left side of most pages displays the navigation icons with a black background.

Home		Ноте	(?) Help	🎗 🛛 Lab Manager User 🗸
Projects	·II¢ Shortcuts	` <u>│</u> ` Latest Notifie	cations	
A Users	\sim	F ⁹ 10/18/18, 8	8:40 AM by LabTech	Ê
ਸਿੰ <u>ਜ</u> ੇ Roles ਜਿੱਸੀ Maintenance	$\left(\begin{array}{c} \mathcal{A}_{+} \end{array} \right) \left(\begin{array}{c} \mathbf{R}_{+} \\ \mathbf{R}_{+} \end{array} \right)$	Image: State	r LabTech could not login because the 8:40 AM by LabTech r LabTech entered an incorrect passwo 8:30 AM by LabTech	ord and wa
System Audit Trail	> Settings > Notifications	> Licenses	8:39 Am by Eublech r LabTech entered an Incorrect passwo 8:39 AM by LabTech r LabTech entered an Incorrect passwo	ord.
	: Information Panel	투 9 10/15/18, 1 Docume	11:26 AM by Scientist ent Status changed. Old Status: In Wor	rk New St
	TOTAL USERS ADMIN USERS	BLOCKED USERS ()	10:46 AM by LabTech ant Status changed. Old Status: In Prog	gress New
		> 1 Users E ^O 10/15/18, 1 Docume	10:26 AM by Scientist ent Status changed. Old Status: In Wor	rk New St
	USED LICENSES ACTIVE PROJECTS	ACTIVE ROLES	9:46 AM by LabTech r LabTech entered an incorrect passwo	ord.
		E 10/15/18, 3	9:44 AM r labted could not login. Invalid uppr	_

User Actions

On the top right, click \checkmark next to your user name:

- Select **Change Date/Time Format** to display the Change Date/Time Format dialog where you can change the format of the date and time display. See Change Date/Time Format on page 25.
- Select **Change Password** to display the Change Password dialog. This option is available for users with the Custom Credentials login mode. See Change Password on page 24.
- Select Logout to log out of the GxP Admin Portal.

Navigation Icons

On the left, the navigation icons display with a black background and provide access to the following workflows.

- Click **Home** to return to the Home page when you navigate away from the Home page.
- Click Projects to display the Project Administration page to manage Projects. Each user with the SoftMax Pro Access permission must be a member of at least one Project. After you add users with the SoftMax Pro Access permission and you define Roles, you assign each user to a Project with a Role. A user can have one Role in each Project and can be a member of multiple Projects. You can assign a user to a different Role in each Project. See Projects on page 42.

- Click **Users** to display the User Administration page to manage users. You grant each user Access permissions. The GxP Admin Access permission grants access to the GxP Admin Portal as an administrator. The System Audit Trail Access permission grants access to the GxP Admin Portal to view the System Audit Trail. The SoftMax Pro Access permission grants access to the SoftMax Pro GxP Software. See Users on page 26.
- Click **Roles** to display the Role Administration page to manage Roles. Use Roles to restrict the responsibilities of the users within the SoftMax Pro GxP Software. Each user with the SoftMax Pro Access permission can be a member of multiple Projects. A user can have one Role in each Project which allows you to assign the same user a different Role with a unique set of permissions in each different Project. See Roles on page 33.
- Click Maintenance to display the Maintenance page. The Maintenance page has four tabs.
 - **Settings**: Use to define auto lock settings, password strength settings, and the administrator contact information. See Maintenance Settings on page 50.
 - **Notifications**: Use to select which events should trigger the notifications that display on the Home page. All events are recorded to the System Audit Trail. See Maintenance Notifications on page 51.
 - Licenses: Use manage the licenses you assign to users with the SoftMax Pro Access permission. See Maintenance Licenses on page 52.
 - User Import: Use to import users from the GxP Admin Software version 2.x into the GxP Admin Software version 3.x database. See Maintenance User Import on page 54.
- Click **System Audit Trail** to display the System Audit Trail page. Use the System Audit Trail to view a list of the events that user actions trigger. When a user has the System Audit Trail Access permission, they can see only the System Audit Trail page. Users with the GxP Admin Access permission automatically receive the System Audit Trail Access permission. See System Audit Trail on page 56.

After you complete the First Steps check list and click Get Started, the Home page displays shortcuts, notifications, and additional information.

Shortcuts

The upper left side of the Home page provides shortcuts to:

- New Users on page 27
- New Roles on page 34
- New Projects on page 43
- Maintenance Settings on page 50
- Maintenance Notifications on page 51
- Maintenance Licenses on page 52

Information Panel

The lower left side of the Home page displays statistics. Each statistic is a link to the workflow where you act on the information.

- Click **Total Users** to display the User Administration page with no filters applied. See Users on page 26.
- Click **Admin Users** to display the User Administration page with filters to display only users with the GxP Admin Access permission.
- Click **Blocked Users** to display the User Administration page with filters to display only users whose software access has been blocked due to too many false login attempts.
- Click **Used Licenses** to display the Maintenance page with the Licenses tab selected where you manage the licenses you assign to users with the SoftMax Pro Access permission. See Maintenance Licenses on page 52.
- Click **Active Projects** to display the Project Administration page with filters to display only active Projects. See Projects on page 42.
- Click **Active Roles** to display the Role Administration page with filters to display only active Roles. See Roles on page 33.

Latest Notifications

The right side of the Home page displays notifications for the events you choose to display. These notifications are similar to your smart phone notifications. When a user triggers an event, the software includes the details in the System Audit Trail.

Use the Notifications tab on the Maintenance page to select which events display as a notification on the Home page. See Maintenance Notifications on page 51.

Note: You cannot change the audit trail. The software permanently records all events in the System Audit Trail even if they do not display as a notification on the Home page.

On the Home page:

- Click on a notification to go to the System Audit Trail page to view additional details.
- Click 🕅 next to a notification to dismiss the notification from the Home page display.
- Click 🗳 on the right to dismiss all notifications from the Home page display.

Change Password

Users with the Custom Credentials login mode use the Change Password dialog to change their password.

Note: Users with the Windows Credentials login mode use the password that is managed via the Windows Active Directory and cannot change their password in the GxP Admin Portal.

Change Password
OLD PASSWORD
\odot
NEW PASSWORD
REPEAT NEW PASSWORD
\odot
Please enter your existing password before setting a new one.
Cancel

To change your password:

- 1. On the top right, click ✓ next to your user name and select **Change Password** to display the Change Password dialog.
- 2. In the **Old Password** field, enter your current password.
- 3. In the **New Password** field, enter your new password.
- 4. In the **Repeat New Password** field, enter your new password a second time to confirm that you enter the password correctly.
- 5. Click Change.

Change Date/Time Format

Use the Change Date/Time Format dialog to define how you want to display the date and time. The database uses the UTC time zone. Your date/time format settings affect the display of the dates on the System Audit Trail page.

Change I	Date,	/Time For	mat	
REGION		FORMAT		UTC TIME
United States (en-US)	~	Short	~	OFF
PREVIEW 10/18/18, 10:28:46 AM				

The Preview area at the bottom of the dialog displays the result of each of the following selections.

To change the date and time format:

- 1. On the top right, click ∨ next to your user name and select **Change Date/Time Format** to display the Change Date/Time Format dialog.
- 2. Click the **Region** drop-down and select a regional date/time format.
- 3. Click the Format drop-down:
 - Select **Short** to display the date and time in the regional format you select.
 - Select **Short + GMT** to display the date and time in the regional format you select plus the offset to the Greenwich Mean Time.
- 4. Click the **UTC Time** to display to have the time display in the regional format you select using the Greenwich Mean Time zone.
- 5. Click Apply.

Chapter 5: Users



E Home			User Administration		? Hel	, R 1	ab Manag	ger User 🗸
Projects	Q Search	Filter					۶,	New User
A Users	USERNAME 👙	FULL NAME	CONTACT INFORM	APPLICATION ACCE	€ STAT	us 😫		
Roles	LabMgr	Lab Manager User	LabMgr@Email.com	Admin, Audit, SMP	Acti	re	Ø	:
& Maintenance	LabTech	Lab Technician User		SMP	Bloc	ked	Ø	:
System Audit Trail	Lab Technician			Admin, Audit	Activ	re	Ø	:
	Scientist	Scientist User	Scientist@Email.com	SMP	Acti	'e	Ø	:

The User Administration page displays the list of users and allows you to manage users.

To manage users:

- 1. Click **Users** on the left to display the User Administration page.
- 2. Click New User to display the New User page. See New Users on page 27.
- 3. Next to a user, click 🖉 to display the Edit User page. See Edit Users on page 30.
- 4. Next to a user, click :

- Select **Reset Password** to reset the user's password. This displays for users with the Custom Credentials login mode. See below.
- Select Deactivate User to make an active user inactive or select Activate User to make an inactive user active.

Note: When you deactivate a user, the software removes all Access permissions and releases the license from a user that had the SoftMax Pro Access permission. You cannot re-activate the user for 30 days.

• Select **User Information** to display user information. See User Information on page 31.

Reset Password

When a user with the Custom Credentials login mode forgets their password, the Reset Password option generates a new alphanumeric string password for the user. The software requires the user to change the system generated password the first time they login.

To reset passwords:

- 1. On the User Administration page, next to a user, click and select **Reset Password**.
- 2. A message displays to confirm your request. You cannot undo this. Click **OK**.
- 3. The User Information dialog displays.
 - Click to copy the user information including the new password to the computer clipboard. Then paste the user information into an email or into a text editor software file and save the file in a location that the user can access.
 - Click Loprint the new user information including the new system generated password.

New Users

Use the New User page to add users.

Select the Login Mode

The first time you click New User, the New User page allows you to select the user's login mode. The next time you click New User, the New User page displays the settings for the login mode you selected for the previous user you added and contains a Change Login Mode button that you use to add a user with the alternate login mode.

ا	Home	New User	(?) Help	R Lab Manager User ∨
Ē	Projects			Save User
8	Users			
Ŕ	Roles	Login Mode		
פצ	Maintenance	& Use Windows Credentials R Use Custom Credentials		
ଔ	System Audit Trail	i Select a login mode for new user.		
		Liss Information		
		oser moniduon		

To select a login mode:

- 1. Click **Users** on the left to display the User Administration page.
- 2. Click **New User** to display the New User page.
 - Click **Use Windows Credentials** to display fields that you use to add a user account that uses your company's Active Directory database for "single sign on" and password management. Windows Credentials login mode is available for the multi computer setup installation only.
 - Click **Use Custom Credentials** to display fields that you use to add a user account that uses the username and password restrictions you define from the GxP Admin Portal.

Windows Credentials Login Mode Users

Windows Credentials login mode users use the username and password settings from your company's Windows Active Directory database.

Note: You must install the SoftMax Pro 7.1.1 GxP Compliance Software Suite using the multi computer setup to use the Windows Credentials login mode. The computer running the GxP Admin Software must have access to the computer that contains your company's Active Directory database.

To add users with the Windows Credentials login mode:

- 1. Click **Users** on the left to display the Users page.
- 2. Click **New User** to display the New User page.
- 3. Click **Use Windows Credentials** to display fields that you use to add a user with the Windows Credentials login mode, if needed.

Note: If you previously added Custom Credential users, click **Change Login Mode** and then click **Use Windows Credentials**.

- 4. In the **Username** field, enter the user's Windows Active Directory username.
- 5. Click **Verify** to link the user's Active Directory account to the GxP Admin Software. The Username field becomes inactive and you cannot edit the Username field.
- 6. The **Full Name** field can display the user's full name from Active Directory or you can enter the full name.
- 7. The **Contact Information** field can display the user's contact information from Active Directory or you can enter the user's contact information. (optional)
- 8. Select Application Access options.
 - Select the **GxP Admin** check box to grant the user Access to the GxP Admin Portal with permission to maintain user information, password strength settings, and Access to the System Audit Trail.
 - Select the **System Audit Trail** check box to grant the user Access to the GxP Admin Portal with permission to view the System Audit Trail.

- 9. Click the **SoftMax Pro Access No Access** → and select a license to grant the user access to the SoftMax Pro GxP Software. Each user with the SoftMax Pro Access permission requires a license and must be a member of a Project. See Maintenance Licenses on page 52 and Projects on page 42.
- 10. Click Save User.

Custom Credentials Login Mode Users

Custom Credentials login mode users use the username and password settings you define from the GxP Admin Portal.

ا	Home	TODO	New	User		? Help Ad Min ∨
Ē	Projects					Save User
R	Users					
Å	Roles		Login Mode (Custom Credentials)		Change Login Mode	
ß	Maintenance		User Information		_	
ଝା	System Audit	t Trail	USERNAME			
			Prease enter a Username A password for the Initial logi	in will be generated automatica	liy.	
			FULL NAME	CONTACT INFORMATION		
			Please enter a Full Name	Please enter Contact Info		
			APPLICATION ACCESS	SOFTMAX PRO ACCESS		
			GxP Admin System Audit Trall	No Access	~	

To add users with the Custom Credentials login mode:

- 1. Click **Users** on the left to display the Users page.
- 2. Click **New User** to display the New User page.
- 3. Click **Use Custom Credentials** to display fields that you use to add a user with the Custom Credentials login mode, if needed.



Note: If you previously added Windows Credential users, click **Change Login Mode** and then click **Use Custom Credentials**.

- 4. In the **Username** field, enter the username.
- 5. In the **Full Name** field, enter the full name.
- 6. In the **Contact Information** field, enter the user's contact information. (optional)
- 7. Select Application Access options.
 - Select the **GxP Admin** check box to grant the user Access to the GxP Admin Portal with permission to maintain user information, password strength settings, and Access to the System Audit Trail.
 - Select the **System Audit Trail** check box to grant the user Access to the GxP Admin Portal with permission to view the System Audit Trail.

- 8. Click the **SoftMax Pro Access No Access** → and select a license to grant the user access to the SoftMax Pro GxP Software. Each user with the SoftMax Pro Access permission requires a license and must be a member of a Project.
- 9. Click Save User.
- 10. The User Information dialog displays. See User Information on page 31.
 - Click I to copy the user information including the new password to the computer clipboard. Then paste the user information into an email or into a text editor software file and save the file in a location that the user can access.
 - Click Loprint the new user information including the new system generated password.

Edit Users

Use the Edit User page to edit users.

ا	Home		Edit Us	er	? Help	8	Lab Manager User 🗸	
Ê	Projects						Save User	
8	Users	Login Mode (Cust	om Credentials)	Change	Login Mode			
Ŕ	Roles	User Information						
פא	Maintenance							
ଟ୍ୟ	System Audit Trail	USERNAME Scientist		CONTACT INFORMATION				
		Scientist User APPLICATION ACCESS GXP Admin System Audit	Trail	Scientist@Email.com SOFTMAX PRO ACCESS SoftMax Pro GXP n.n The license stoys with the user for 23	✓ days and			
				cannot be changed within this time p	eriod.			

To edit users:

- 1. Click **Users** on the left to display the Users page.
- 2. Next to a user, click 🖉 to display the Edit User page.
- To change a Custom Credentials login mode user to a Windows Credentials login mode user or to change a Windows Credentials login mode user to a Custom Credentials login mode user, click Change Login Mode and then click either Use Windows Credentials or click Use Custom Credentials.
- 4. In the **Username** field, enter the username. You cannot change the username for a Windows Credentials login mode user.
- 5. In the **Full Name** field, enter the full name.
- 6. In the **Contact Information** field enter the user's contact information.

- 7. Select Application Access options.
 - Select the **GxP Admin** check box to grant the user Access to the GxP Admin Portal with permission to maintain user information, password strength settings, and Access to the System Audit Trail.
 - Select the **System Audit Trail** check box to grant the user Access to the GxP Admin Portal with permission to view the System Audit Trail.
- 8. Click the **SoftMax Pro Access No Access** → and select a license to grant the user access to the SoftMax Pro GxP Software. Each user with the SoftMax Pro Access permission requires a license and must be a member of a Project. See Maintenance Licenses on page 52 and Projects on page 42.
- 9. Click Save User.

User Information

The User Information dialog displays user information.

User Informatio	on
Lab Manager 2 Laboratory Manager	
CONTACT INFORMATION LabManager@Email.com LOGIN MODE Custom Credentials TEMPORARY PASSWORD d457c5590fbc4c97ac56 (i) STATUS Active APPLICATION ACCESS Admin, Audit, SMP	
Close	

To view user information:

- 1. Click **Users** on the left to display the User Administration page.
- 2. Next to a user, click and select **User Information** to display the User Information dialog.
- 3. Click to copy the user information to the computer clipboard from where you can paste the information into an email or text file.
- 4. Click 💾 to print the user information.

Chapter 6: Roles

The Role Administration page displays the list of Roles. Use the Role Administration page to manage Roles. Use Roles to define the list of permissions that grant a user access to the workflows in the SoftMax Pro GxP Software. You assign a role to a user when you manage Projects from the Project Administration page. See Projects on page 42.

Each user with the SoftMax Pro Access permission can be a member of multiple Projects. The user has one Role in each Project. This allows you to assign the same user a different Role with a unique set of permissions in each Project.

The software installation includes three default Roles. You can edit the default Roles or copy the default Roles to create your own similar Roles. See Default Roles on page 63.

Home TODO	Role Administration					Jser2 ∨
Projects		Filter				ew Role
A Users	ROLE NAME		PERMISSIONS	PROJECTS 👙	STATUS 😝	
Roles	Lab Manager	Standard Workflow	26	0	Active 🖉	:
B Maintenance	Lab Technician	Standard Workflow	12	0	Active 🖉	:
System Audit Trail	Scientist	Standard Workflow	42	0	Active 🖉	:

To manage Roles:

- 1. Click **Roles** on the left to display the Role Administration page.
- 2. Click New Role to add a new Role. See New Roles on page 34.
- 3. Next to a Role, click or to display the Edit Role page. See Edit Roles on page 35.
- 4. Next to a Role, click 🔹 :
 - Select **Copy Role** to display the Copy Role page. See Copy Roles on page 36.
 - Select **Deactivate Role** to make an active Role inactive or **Activate Role** to make an inactive Role active.
 - Select **Assigned Permissions** to display the Assigned Permission dialog that displays the permission that are granted to users with the Role. See Assigned Permissions on page 37.
 - Select **Assigned Projects** to display the Assigned Projects dialog that displays the list of Projects that include users with the Role. See Project Information on page 48.

New Roles

? Help ♀ Lab Manager User ∨ 🔓 Home New Role Projects Role Name 🖉 ROLE NAME A Users Access Select All Roles DESCRIPTION nported Open In Work Open Review Pending Ŗ Mainte Open Released Open Outdated eviewed System Audit Trail Open Approval Pending Open Approved Progress anceled ок Select All Workflow O Set Review Pending O Release Protocol Set Reviewed Generate compliance data Set Outdated Set Approval Pending Set Approved Set Canceled Document Editing Select All Read empty plates/cuvettes O Edit reader settings Edit notes text Overwrite plate/cuvette data Edit graphs, summaries and reductions O Function editor

Use the New Role dialog to add Roles.

To add Roles:

- 1. Click **Roles** on the left to display the Role Administration page.
- 2. Click **New Role** to display the New Role page.
- 3. In the **Role Name** field, enter the Role name.
- 4. In the **Description** field, enter the Role description and click **OK**.
- 5. Select the check box for each permission to grant to the users you assign to the Role. See Role Permissions on page 39.
- 6. Click Save Role.

Edit Roles

Home			Edit Role		? Help ♀ U2 Lab Manager ∨
Projects	Scientist 🖉				Save Role
A Users				0	
Roles		Document Access		Select A	11
Maintenance		Open Imported	Open In Work	Open Review Pending	
		Open Reviewed	Open Released	Open Outdated	
System Audit Trail		Open In Progress	Open Approval Pending	Open Approved	
		Open Canceled			
		Document Workflow		Select A	11
		Set Review Pending	Set Reviewed	Release Protocol	
		Set Outdated	Generate compliance data	Set Approval Pending	
		Set Approved	Set Canceled		
		Document Editing		Select A	n
		Edit reader settings	Rdlt notes text	👩 Read empty	

Use the Edit Role page to edit Roles.

To edit Roles:

- 1. Click **Roles** on the left to display the Role Administration page.
- 2. Next to a Role, click 🖉 to display the Edit Role page.
- 3. On the top left, click of to edit the **Role Name** and **Description**.
- 4. Select the check box for each permission to grant to the users you assign to the Role. See Role Permissions on page 39.
- 5. Click Save Role.

Copy Roles

Use the Copy Role page to create copies of existing Roles to expedite the creation of similar Roles.



To copy Roles:

- 1. Click Roles on the left to display the Role Administration page.
- 2. Next to a Role, click and select **Copy Role** to display the Copy Role page.
- 3. In the **Role Name** field, enter the Role name.
- 4. In the **Description** field, enter the Role description and click **OK**.
- 5. Select the check box for each permission to grant to the users you assign to the Role. See Role Permissions on page 39.
- 6. Click Save Role.
Assigned Permissions

	S
Scientist	
43 PERMISSIONS	
Add folder	
Add/Delete groups	
Add/Modifly statements	
Assign plate layouts	
Change Auto Export settings	
Change Auto Save settings	
Create/Save as data document	
Create/Save as protocol	
Delete document	

The Assigned Permissions dialog displays the permissions you assign to a Role.

To view assigned permissions:

- 1. Click **Roles** on the left to display the Role Administration page.
- 2. Next to a Role, click and select **Assigned Permissions** to display the Assigned Permissions dialog.
- 3. Click to copy the permissions to the computer clipboard from where you can paste the information into an email or text file.
- 4. Click 📕 to print the permissions.

Assigned Projects

The Assigned Projects dialog displays the list of Projects that use the Role and the number of users in the Role for each Project.

Assigned Pi	rojects
Scientist	
2 ASSIGNED PROJECTS	
Inactive Project	1 Users
Workflow Project	1 Users
Close	

To view assigned Projects:

- 1. Click **Roles** on the left to display the Role Administration page.
- 2. Next to a Role, click and select **Assigned Projects** to display the Assigned Projects dialog.
- 3. Click copy the assigned Projects list to the computer clipboard from where you can paste the information into an email or text file.
- 4. Click 💾 to print the assigned Projects list.

Role Permissions

In the GxP Admin Portal, you assign permissions to a Role. When you assign users to a Project, you select each user's Role in the Project. For each Project, the user can have one Role that permits the user to perform their required tasks, and no more. You can assign each user to multiple Projects and the user can have a different Role in each Project. This provides extensive and highly granular administrative control over the user's activities in the SoftMax Pro GxP Software.

Role permissions are grouped into the following categories:

Document Access Permissions

Use Document Access permissions to manage what users can do with documents.

- **Open Imported** Allows users to open documents that have the status: Imported.
- **Open In Work** Allows users to open documents that have the status: In Work.
- **Open Review Pending** Allows users to open documents that have the status: Review Pending.
- **Open Reviewed** Allows users to open documents that have the status: Reviewed.
- **Open Released** Allows users to open documents that have the status: Released.
- **Open Outdated** Allows users to open documents that have the status: Outdated.
- **Open In Progress** Allows users to open documents that have the status: In Progress.
- **Open Approval Pending** Allows users to open documents that have the status: Approval Pending.
- **Open Approved** Allows users to open documents that have the status: Approved.
- **Open Canceled** Allows users to open documents that have the status: Canceled.

Document Workflow Permissions

Use Document Workflow permissions to manage what users can do in the document workflow.

- Set Review Pending Allows users to set documents to the status: Review Pending.
- Set Reviewed Allows users to set documents to the status: Reviewed.
- **Released Protocol** Allows users to release a document as a protocol which sets the document status to Released.
- Set Outdated Allows users to set documents to the status: Outdated.
- Generate Compliance Data Allows users to generate compliance data. For example, (using the system default Roles) this user would be the Lab Technician who can open a protocol document with the status Released. The software prompts the user to save the document as a data document and assigns the document the status In Progress. The user then runs the assay to generate compliance data.
- Set Approval Pending Allows users to set documents to the status: Approval Pending.
- Set Approved Allows users to set documents to the status: Approved.
- Set Canceled Allows users to set documents to the status: Canceled.

Document Editing Permissions

Use Document Editing permissions to manage what users can do with documents.

- Edit Reader Settings Allows users to define microplate reader acquisition settings.
- Edit Notes Text Allows users to edit Note sections in the experiments in documents.
- Read Empty Plates/Cuvettes Allows users to read empty Plate sections and empty Cuvette Set sections.
- Function Editor Allows users to create custom curve fit functions.
- **Overwrite Plate/Cuvette Data** Allows users to overwrite the data in Plate sections and Cuvette Set sections.
- Edit Graphs, Summaries, and Reductions Allows users to edit Graph sections and to edit Summary formulas and Reduction formulas.
- Assign Plate Layouts Allows users to assign plate layouts.
- Edit Sample and Group Information Allows users to edit sample and Group section information.
- Add/Delete Groups Allows users to add and delete Group sections from the experiments in documents.
- Edit Formulas Allows users to edit formulas in documents.
- Edit Print Options Allows users to edit the print options.
- Lock/Unlock Sections Allows users to lock and unlock sections in the experiments in documents.
- Mask/Unmask Wells Allows users to mask and unmask wells.
- Create/Save As Data Document Allows users to save data documents to the database.
- Create/Save As Protocol Allows users to save protocol documents to the database.
- Change Auto Save Settings Allows users to change Auto Save settings.
- Change Auto Export Settings Allows users to change Auto Export settings.

Statements & Signatures Permissions

Use Statement and Signatures permissions to manage what users can do with the statements.

- Add/Modify Statements Allows users to edit statements.
- Sign Statements Allows users to sign statements.
- Revoke Own Signature Allows users to remove only their own signatures from statements.
- Revoke Any Signature Allows users to remove any signature from statements.

Document Management Permissions

Use Document Management permissions to manage what users can do with the documents in the database.

- Rename Document Allows users to rename documents.
- Move Document Allows users to move documents to different folders in the database.
- Delete Document Allows users to delete documents from the database.
- Unlock Document Allows users to unlock documents.
- Import Documents Allows users to import documents into the database.
- **Export Documents** Allows users to export documents out of the database.

Folder Management Permissions

Use Folder Access permissions to manage what users can do with the document storage folders in the database.

- Add Folder Allows users to create new folders in the database.
- Rename Folder Allows users to rename folders in the database.
- Move Folder Allows users to move folders in the database.
- Delete Empty Folder Allows users to delete empty folders from the database.
- **Hide/Unhide Folder** Allows users to hide and unhide folders and to use hidden folders in the database.

Instrument Permissions

Use Instrument permissions to manage what users can do with the microplate reader.

- Lock/Unlock Instrument Allows users of the SpectraMax[®] iD3 Multi-Mode Microplate Reader and the SpectraMax[®] iD5 Multi-Mode Microplate Reader to lock and unlock the instrument touchscreen.
- Instrument Simulator Allows users to simulate the connection between the SoftMax Pro GxP Software and an instrument.

Chapter 7: Projects

шь



The Project Administration page displays the list of Projects. Use the Project Administration page to manage Projects. Each user with the SoftMax Pro Access permission must be a member of at least one Project before they can login to the SoftMax Pro GxP Software. Each user with the SoftMax Pro access permission can be a member of multiple Projects and the user can have a different Role in each Project.

	Note: P	Prerequisites: Yo	ou must define u	users and R	oles before yo	u create Project	5.
۵	Home			Project Adminis	stration	? Help A Lab Manag	er User 🗸
Ê		Q Search	Filter			Net Ne	w Project
8	Users	PROJECT NAME	DESCRIPTION	USERS 🖨	ROLES 👙	STATUS 😝	
Ŕ	Roles	Valldation Project	Validation Protocol Te	2	2	Active 🖉	:
Ŗ	Maintenance	Workflow Project	Document Workflow	3	3	Active 🖉	:
ଝା	System Audit Trail						

To manage Projects:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Click **New Project** to display the New Projects page. See New Projects on page 43.
- 3. Next to a Project, click or to display the Edit Project page. See Edit Projects on page 44.
- 4. Next to a Project, click :
 - Select **Copy Project** to display the Copy Project page. See Copy Projects on page 45.
 - Select **Deactivate Project** to make an active Project inactive or select **Activate Project** to make an inactive Project active.
 - Select **Project Information** to display the Project Information dialog. See Project Information on page 48.

New Projects

Use the New Project page to add Projects.

ຝ	Home	New Project	? Help	8 Lab Manager User ∨
Ê	Projects	Project Name 🖉		Save Project
Я	Users		0 Add1	leare
Ŕ	Roles		/ /+ Aud (
פא	Maintenance	Please enter an optional description No users added yet.		
ଷ୍	System Audit Trail			
		ок		

To add Projects:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Click **New Project** to display the New Project page.
- 3. In the **Project Name** field, enter the Project name.
- 4. In the **Description** field, enter the Project description and then click **OK**.
- 5. Click **Add Users** to display the Assign Users to Project Role dialog where you add users to the Project and assign users to a Role in the Project. See Assign Users to Project Roles on page 46.
- 6. After you add users, click **Save Project** to save the project.

Edit Projects

- Home			Edit Project		? Help & La	ib Manager User 🗸 🗸
Projects	Workflow Project	Ŷ				Save Project
A Users						
Roles						
Maintenance	_	USERNAME	ROLE 😫	USER STAT	\$	
		LabMgr	Lab manager	Active	:	
System Audit Trail		LabTech	Lab technician	Blocked	:	
		Scientist	Scientist	Active	:	
			3 users in total.			

Use the Edit Project page to edit Projects.

To edit Projects:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Next to a Project, click 🖉 to display the Edit Project page.
 - On the top left, click Ø to edit the **Project Name** and **Description**.
 - Click **Add Users** to display the Assign Users to Project Role dialog where you add users to the Project and assign users to a Role in the Project. See Assign Users to Project Roles on page 46.
 - Click next to the Add User button to display a check box next to each username. Select the check box for each user to remove from the Project and then click to remove the users you select from the Project.
- 3. Next to a username, click :
 - Select **Change Role** to display the Change Role dialog where you change the user's Role in the Project. See Change Role on page 47.
 - Select **Remove User** to remove the user from the Project.
- 4. Click Save Project.

Copy Projects

Home		Copy Project	⑦ Help
Projects	Workflow Project Copy 🔗 Document Workflow		Save Project
A Users	PROJECT NAME Workflow Project Copy		₩ Ø A. Add Users
हिं <u>ः</u> Roles	Description Document Workflow	ROLE 🔶	USER STAT 🗧
॰ Q System Audit Trail		Lab manager	Active
		Lab technician	Blocked
	ОК	Scientist	Active
		З users in total.	

Use the Copy Project page to copy Projects.

To copy Projects:

- 1. Click **Projects** on the left to display the Projects page.
- 2. Next to a Project, click and select **Copy Project** to display the Copy Project page.
- 3. In the **Project Name** field, enter the Project name.
- 4. In the **Description** field, enter the Project description and click **OK**.
- 5. Click **Add Users** to display the Assign Users to Project Role page where you add users to the Project and assign users to a Role in the Project. See Assign Users to Project Roles on page 46.
- 6. Next to a username, click :
 - Select **Change Role** to display the Change Role dialog where you change the user's Role in the Project. Change Role on page 47.
 - Select **Remove User** to remove the user from the Project.
- Click [→] P next to the Add User button to display a check box next to each username.
 Select the check box for each user to remove from the Project and then click [□] to remove the users you select from the Project.
- 8. Click Save Project.

Assign Users to Project Roles

_

Use the Assign Users to Project Role dialog to add users to the Project and to assign users to a Role in the Project. You must define Users and Roles before you can assign users to a Project.

	Assign Use	ers to Project Role
SELECT FROM ALL USERS		SELECTED USERS
Q Search		LabMgr
LabMgr	\bigcirc	
LabTech	\oplus	
Lab Technician	\oplus	
Scientist	\oplus	
SELECT FROM ALL ROLES		
Please select a Role	× 0	Choose a Role for the selected users.

To add users to a Project and assign users a Role in the Project:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Click **New Project** to display the Add Projects page or click next to a Project to display the Edit Projects page. See New Projects on page 43 and Edit Projects on page 44.
- 3. Click Add Users to display the Assign Users to Project Role dialog.
- 4. In the **Select From All Users List**, next to each user to assign to the first Role, click to each user to move the name to the Selected Users column.
- 5. Click the **Please Select A Role V** and select a Role for the users.
- 6. Click Assign.
- 7. Click **Add Users** and repeat steps 4-6 to add additional users to a second Role in the Project, and so on for additional Roles, as needed.

Change Role

Use the Change Role dialog to change the user's Role in a Project.

	Char	ige Rol	le	
USERNAME User2 FULL NAME U2 Lab Manager CURRENT PROJECT ROLE Lab Manager				
SELECT FROM ALL ROLES Please select a Role		~		
	Cancel		ОК	

To change user Role assignments:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Click **New Project** to display the Add Projects page or click *Project* to a Project to display the Edit Project page. See New Projects on page 43 and Edit Projects on page 44.
- 3. Next to a username, click and select **Change Role** to display the Change Role dialog.
- 4. Click the **Please Select A Role ∨** and select a Role for the user.
- 5. Click OK.

Project Information

Proje	ect Information
Workflow Project Document Workflow 1 USERS ASSIGNED	
LabMgr	Lab manager
	Close

The Project Information dialog displays the list of users and their Role assignment in the Project.

To display assigned projects:

- 1. Click **Projects** on the left to display the Project Administration page.
- 2. Next to a Project, click and select **Project Information** to display the Project Information dialog.
- 3. Click to copy the information to the computer clipboard from where you can paste the information into an email or a text file.
- 4. Click 💾 to print the information.

Chapter 8: Maintenance



Use the Maintenance page to define password strength settings, to define which events display a notification on the Home page, to manage the licenses you assign to users with the SoftMax Pro Access permission, and to import legacy usernames.

۵	Home	Maintenance	? Help ႙ Lab Manager User ∨
Ē	Projects	SETTINGS NOTIFICATIONS LICENSES USER IMPORT	
8	Users		Save
Ŕ	Roles	General Settings	
		Auto Lock An inactive user will be locked after this number of minutes (1 - 60). This setting is SMP only.	ON • 5 ¢
ଗ୍ର	System Audit Trail	Jenny Administrator 867-5309	
		Password Settings ①	
		Minimum Password Length From minimum 6 to maximum 30 characters.	6 A _V
		Requires Uppercase Letters	
		Requires Lowercase Letters	
		Reautres Diatis	ON O

The Maintenance page has four tabs.

- Select the **Settings** tab to enter administrator contact information, define auto lock settings, and define password strength options. See below.
- Select the **Notifications** tab to define which events should trigger a notification to appear on the Home page. See Maintenance Notifications on page 51.
- Select the **Licenses** tab to manage the licenses you assign to users with the SoftMax Pro Access permission. See Maintenance Licenses on page 52.
- Select the **User Import** tab to import GxP Admin Software version 2.x usernames into the GxP Admin Software version 3.x database. See Maintenance User Import on page 54.
 - **Note:** Windows Credentials login mode users use your company's Active Directory password settings.

Maintenance Settings

Use the Settings tab on the Maintenance page to define auto lock settings for the SoftMax Pro GxP Software, administrator contact information, and password strength options. The default password strength settings require Custom Credentials login mode user passwords to contain at least 10 characters, both uppercase and lowercase letters, at least one digit, and at least one special character.

To manage maintenance settings:

- 1. Click **Maintenance** on the left to display the Maintenance page with the **Settings** tab selected by default.
- 2. Click the **Auto Lock (e**) to display **(*•)**. Enter a duration between 1 and 60 minutes to disable a user's session if they are inactive in the SoftMax Pro GxP Software for the duration you set. This does not affect the GxP Admin Portal.
- 3. In the **Administrator Contact Information** field, enter the name and contact information for the person who can reset a user's password and can unblock a user account when the user enters too many failed login attempts.
- 4. In the Password Settings section, set each of the following to to activate the feature:
 - Use the **Minimum Password Length** field to define the minimum number of characters for user passwords. Enter between 6 and 30 characters.
 - Use the **Requires Uppercase Letters** setting to require users to include upper case letters in their password.
 - Use the **Requires Lowercase Letters** setting to require users to include lower case letters in their password.
 - Use the **Requires Digits** setting to require users to include numbers in their password.
 - Use the **Special Characters** setting to require users to include special characters in their password.
 - Use the **Password Has To Differ From Username** setting to prevent users from using their username as their password.
 - Use the **Password Aging** setting to force users to change their password every x number of days. Enter the number of days that can elapse before the user must change their password, between 30 and 365 days.
 - Use the **Password History** setting to prevent users from reusing their previous x number of passwords. Enter the number of passwords that cannot be repeated, between 2 and 20.
 - Use the Max. Number of False Login Attempts setting to have the system block a user account after a number of incorrect password entries. Enter the number of allowable attempts between 3 and 10.

Maintenance Notifications

Use the Notifications tab on the Maintenance page to select which events display a notification on the Home page each time a user triggers the event.

ا	Home	TODO			Maintenance	? Help	🎗 🛛 Lab Manager User 🗸
ê	Projects		SETTING	S NOTIFICATIONS LICENSES	USER IMPORT		
R	Users						Save
Å	Roles		Notif	flable Events (j			
			•	Document - Released Protocol Created	User created a Released protocol from a Reviewed document.		
ସ୍ତ୍ର	System Aud	it Trail	Ø	Document - Status Changed	User changed the status of a document.		
			0	Project - Edited	User changed a Project.		
			0	Role - Edited	User changed a Role.		
			0	System - Server License Installed	User added additional SoftMax Pro Software access licenses.		
			0	System - Settings Changed	User changed settings on the GxP Admin Portal.		

To select the events to display notifications:

- 1. On the Maintenance page, select the **Notifications** tab.
- 2. Select the check box for each event that you want to display a notification on the Home page.
- 3. Click Save.

Maintenance Licenses

Use the Licenses tab on the Maintenance page to activate the licenses for users with the SoftMax Pro Access permission.

Each user to whom you grant the SoftMax Pro Access permission requires a license.

🔓 Home				Maintenance		? Help	R Lab Manager User ∨
Projects	SETTINGS	NOTIFICATIONS	LICENSES	USER IMPORT			
R Users			_				
Roles	License Co	de (i)					
🔑 Maintenance	Enter Licer		Get Licenses				
System Audit Trail							
	Application	Access Ve	rsion	Total Licenses	Used Licenses	Available	Licenses
	Application SoftMax Pr	o GXP 7.1	rsion	Total Licenses	Used Licenses	Available 20	Licenses
	Application SoftMax Pr	o GxP 7.1	rsion	Total Licenses	Used Licenses	Available 20	Licenses
	Application SoftMax Pr	o GXP 71	rsion	Total Licenses	Used Licenses	Available 20	Licenses
	Application SoftMax Pr	ı Access Ve o GXP 7.1	rsion	Total Licenses	Used Licenses	Available 20	Licenses
	Application SoftMax Pr	I Access Ve	rsion	Total Licenses	Used Licenses	Available 20	Licenses

When you use upgrade licenses for multiple software versions of the GxP Admin Software and you end up with more users than licenses, use the License Upgrade dialog to assign applicable licenses to users. See License Upgrade on page 53.

Computer Has Internet Access

To activate licenses on a computer with Internet Access:

- 1. On the Maintenance page, select the Licenses tab.
- 2. In the **License Code** field, enter the license code.
- 3. Click Get Licenses.

Computer Does Not Have Internet Access

If the computer does not have Internet access:

- 1. In the **License Code** field, enter the license code.
- 2. Click Get Licenses.
- 3. Click **Export** and save the file that the system creates to a flash drive.
- 4. Take the flash drive to a computer that has Internet Access and go to https://smplicensing.moleculardevices.com/gxpreg.
- 5. Follow the instructions to activate the license and save the Import file to the flash drive.
- 6. Take the flash drive back to the computer running the GxP Admin Portal and click **Import**.

License Upgrade

When you use upgrade licenses for multiple versions of the GxP Admin Software and you end up with more users than available licenses, use the Licenses Upgrade dialog to resolve license issues. Each user with the SoftMax Pro Access permission requires a license.

License Upgrade	
License Conflict ① SoftMax Pro GxP 15 users assigned, 14 licenses in total.	
Q Search	
LabMgr	SoftMax Pro GxP
LabTech	SoftMax Pro GxP
Scientist	SoftMax Pro GxP V
Rebecca	SoftMax Pro GxP · V
Alicia	SoftMax Pro GxP V
Ted	SoftMax Pro GxP V
Cancel Assig	n

To resolve license issues:

- 1. Next to each user, click the **SoftMax Pro GxP (n.n)** V and select a license.
- 2. Click Assign.

Maintenance User Import

Use the User Import tab on the Maintenance page to import the names of the users you created in GxP Admin Software version 2.x into the GxP Admin Software version 3.x database. Use the GxP Admin EDB Converter Tool to convert the version 2.x *.edb files to the version 3.x *.edbexp files before you import usernames into the 3.x database. See Converting Legacy User Files on page 16.

Ē

Note: This process imports usernames, passwords, and the GxP Admin Access permission, when applicable.

After you complete the import, you must assign each user with the SoftMax Pro Access permission a license and to at least one Project.



You repeat the following steps for each .edbexp file you create with the GxP Admin EDB Converter Tool.

To import users:

- 1. On the Maintenance page, select the **User Import** tab.
- 2. Click Import From File to display the Open dialog.
- 3. Navigate the folder hierarchy and select an .edbexp file to import.
- 4. Click Open to display the Import Legacy Users dialog.

Select Users to Import

Use the Import Legacy Users dialog to select the individual usernames to import into the GxP Admin Software version 3.x database.

	Import Legacy Users
SELECT LEGACY USERS	SELECTED USERS
Q Search	
Select All	\oplus
Rebecca	\oplus
Alicia	\oplus
	Cancel Import

To select the users to import:

- 1. In the Select Legacy Users column, next to each user to import, click \oplus to move the name to the Selected Users column.
- 2. Click Import.

Chapter 9: System Audit Trail

Use the System Audit Trail page to view the events that user actions trigger. You cannot edit or delete System Audit Trail entries. You can change the display of the page and you can filter the list of events to hide events from the display. The database uses the UTC time zone. Your date/time format settings affect the display of the dates on the System Audit Trail page. See Change Date/Time Format on page 25.

<	< Home			System Audit Trail		? Help	8	Lab Manager User 🗸
	Layout				C Refresh	Add Note		Create Report
	DATE/TIME	EVENT 😫	USERNAME	DESCRIPTION	DOCUMENT			√ Filter Settings
	11/1/18, 1:24 PM	User - Logged In	LabMgr	The User LabMgr has logged in.				No filters applied.
	11/1/18, 1:16 PM	User - Logged Out	LabMgr	The User LabMgr has logged out.				
	11/1/18, 12:26 PM	User - Logged In	LabMgr	The User LabMgr has logged in.				
	11/1/18, 12:11 PM	User - Logged Out	LabMgr	The User LabMgr has logged out.				
	11/1/18, 11:41 AM	Project - Added	LabMgr	User added a Project. Project Name: Wor				
	11/1/18, 11:28 AM	User - Logged In	LabMgr	The User LabMgr has logged in.				
	10/31/18, 3:39 PM	User - Logged Out	LabMgr	The User LabMgr has logged out.				
	10/31/18, 3:24 PM	User - Logged In	LabMgr	The User LabMgr has logged in.				
	10/31/18, 3:07 PM	User - Logged Out	LabMgr	The User LabMgr has logged out.				
	10/31/18, 2:34 PM	User - Logged In	LabMgr	The User LabMgr has logged in.				

Click **System Audit Trail** on the left to display the System Audit Trail page.

- Click **Layout** to display the Custom Layout dialog where you define which columns of data display. See System Audit Trail Layout on page 57.
- Click **Refresh** to update the display.
- Click Add Note to display the Add Note To Audit Trail dialog where you add a note. See System Audit Trail Notes on page 59.
- Click **Create Report** to display the Create Report dialog where you create a report. See System Audit Trail Report on page 60.
- Click **Filter Settings** to filter the list of events that display. See System Audit Trail Filters on page 58.
- Click **Home** to return to the Home page. See Home on page 21.

System Audit Trail Layout

Use the Custom Layout dialog to define which columns of data display in the System Audit Trail page.



The System Audit Trail can display six columns of data. The Data/Time column must remain visible and must be located in the far left column.

- 1. Click **System Audit Trail** on the left to display the System Audit Trail page.
- 2. Click **Layout** to display the Custom Layout dialog.
 - The **Visible** area on the left displays the data columns that display in the System Audit Trail page. Each visible column has a light background and appears in the sequence in which they display.
 - The **Hidden** area on the right displays the data columns that do not display in the System Audit Trail page. Each hidden column has a shaded background.
- 3. Click on a column to display a blue border around the column, then click **Move Left** or click **Move Right** to change the column sequence and to move columns to the Visible area or to the Hidden area. Columns in the Visible area display in the System Audit Trail page in the sequence that displays in the Custom Layout dialog.
- 4. Click **OK** to display the System Audit Trail page.

System Audit Trail Filters

<	Home			System Audit Trail	R Lab Manager User ∨
	Layout			C Refresh & Add Note	Create Report
	DATE/TIME	EVENT 👙	USERNAME 👙		USER FILTER (0/5)
	10/18/18, 2:52 PM	EDB-File imported	LabMgr	Q Filter.	
	10/18/18, 2:52 PM	Admin permission	LabMgr		
	10/18/18, 2:52 PM	User created	LabMgr	C Lab Manager 2	EVENT FILTER (0/5)
	10/18/18, 2:52 PM	EDB-File imported	LabMgr	O LabMgr	DOCUMENT FILTER (0/5) DATE FILTER (0/1)
	10/18/18, 2:34 PM	Login	LabMgr	O Rebecca	
	10/18/18, 2:29 PM	Logout	LabMgr	O Scientist	
	10/18/18, 12:59 PM	License assigned	LabMgr		
	10/18/18, 12:59 PM	Admin permission	LabMgr		
	10/18/18, 12:59 PM	User created	LabMgr	Filters Applied: 1109 events will display.	
	10/18/18, 12:40 PM	Login	LabMgr	Remove Filter	Close

Use the Filter Settings dialog to limit the events that display in the System Audit Trail page. The Filter Settings dialog has four tabs and you can set up to five filters per tab.

To filter the list of events that display on the System Audit Trail page:

- 1. Click **System Audit Trail** on the left to display the System Audit Trail page.
- 2. Click **Filter Settings** to display the Filter Settings dialog.
- 3. The Filter Settings dialog has four tabs. After you select the tab, select up to five check boxes on the left. The fields on the right display the list of filters you select.
 - Use the **User Name** tab to select up to five users to limit the list of events that display on the System Audit Trail page to only events triggered by the users you select.
 - Use the **Events** tab to select up to five events to limit the list of events that display on the System Audit Trail page to only the events you select. See Audit Events on page 66.
 - Use the **Documents** tab to select up to five documents to limit the list of events that display on the System Audit Trail page to only the events related to the documents you select.
 - Use the **Date** tab to select up to five dates to limit the list of events that display on the System Audit Trail page to only events that occurred on the dates you select.
- 4. Click **Apply** to display the System Audit Trail page with the list of events filtered to the display only the events that meet your filter criteria.
- 5. Click **Remove Filter** to remove all filters.

System Audit Trail Notes

Use the Add Note To System Audit Trail dialog to add a note to the System Audit Trail.

Add Note to System Audit Trail
ENTER TEXT FOR NOTE
• The note is added to the System Audit Trail and becomes part of the permanent record.
Cancel Add

To add a note to the System Audit Trail:

- 1. On the left, click **System Audit Trail** to display the System Audit Trail page.
- 2. Click Add Note to display the Add Note To System Audit Trail dialog.
- 3. In the Enter Text For Note field, enter the note text.
- 4. Click **Add** to add the note to the audit trail.

System Audit Trail Report

Use the Create Report dialog to create a System Audit Trail report file that you can save with a .pdf file format or a .csv file format.

Report Settings		
FILE FORMAT		
PDF	~	
PAPER FORMAT		
Α4	~	
LAYOUT OPTIONS		
Archive Layout		
O Binding Margin		

To create a System Audit Trail report:

- 1. On the left, click **System Audit Trail** to display the System Audit Trail page.
- 2. Click **Create Report** to display the Create Report dialog.
- 3. To create a report with the .pdf file extension, click the **File Format** V and select **PDF**:
 - a. Click the **Paper Format** \checkmark and select the paper format.
 - b. Select the **Archive Layout** check box to create a report in a list format. Leave this check box clear to create a report in a table format.
 - c. Select the **Binding Margin** check box to justify the text to accommodate a binding. Leave this check box clear to center the report on the page.
- 4. To create a report with the .csv file extension, click the **File Format** and select **CSV**. The **Delimiter** field displays **CSV**.
- 5. Click **Generate**, the report generates and the Open dialog displays.
- 6. Navigate to where you want to save the report file and click **Save**.



Chapter 10: Logging In To The SoftMax Pro Software - GxP Edition

When you start the SoftMax Pro Software - GxP edition, the SoftMax Pro GxP Log On dialog displays.

🚺 SoftMax Pro GxP – Log On			×
STICFR Part	<user id=""></user>		
CARP OF CONTRACT O	Show Password		
	Contact:		
DEVICES			
		Next	xit SMP

To log on to the SoftMax Pro Software - GxP edition:

- 1. In the **<User ID>** field, enter your user name.
- 2. In the **<Password>** field, enter your password (case sensitive).
- 3. Click **Next** to display the Please Select a Project dialog.

🔽 SoftMax Pro GxP – Log C	Dn Please select a project:	×
	Workflow Project	
SoftMax Pro GXP ONPLANE	Validation Project Role: Lab manager	
DEVICES		
	Log In	Back

4. Select a Project and click Log In.

Logging On From Within The Software

In the Ribbon, on the GxP tab, use the icons to log out of the software and then log on to a different Project or as a different user.

- 1. In the Ribbon, select the **GxP** tab.
- 2. Click ^O User Log Off to log off. The SoftMax Pro GxP Log On dialog displays.
- 3. Follow the steps above to log on.

SpectraMax iD3 or iD5 Considerations

In the Ribbon, on the GxP tab, users with appropriate permission can use the following icons to lock and unlock the instrument touchscreen:

- Click **GxP Mode On** to lock the instrument touchscreen and operate the instrument from the computer running the SoftMax Pro Software in GxP mode. This locks the instrument touchscreen for all users and you must operate the instrument from a computer running the SoftMax Pro Software GxP edition.
- Click **GxP Mode Off** to release the lock from the instrument touchscreen and allow users to use the instrument touchscreen to run experiments.

Note: The instrument remains locked until the user with the appropriate permission

clicks **GxP Mode Off** to stop the GxP mode. You cannot use the Instrument Connection dialog to disconnect from a SpectraMax iD3 and SpectraMax iD5 that is locked in GxP mode.

Appendix A: Default Roles



Use the GxP Admin Software default Roles to assign users permissions to perform the tasks that correspond to each document status. You can create your own Roles and you can edit the default Roles.

Default	Role	Permi	ssions
Derduit	NOIC	I CITILI	3310113

Permission Group	Permission Name	Lab Mgr	Lab Tech	Scientist
Document Access	Open Imported			\checkmark
	Open In Work			\checkmark
	Open Review Pending	~		
	Open Reviewed	~		\checkmark
	Open Released	~	~	\checkmark
	Open Outdated	~		\checkmark
	Open In Progress		~	
	Open Approval Pending	~		\checkmark
	Open Approved	~		\checkmark
	Open Canceled	~		\checkmark
Document Workflow	Set Review Pending			\checkmark
	Set Reviewed	~		
	Release Protocol	~		
	Set Outdated	~		
	Generate Compliance Data		~	
	Set Approval Pending		~	
	Set Approved	~		
	Set Canceled	\checkmark	\checkmark	\checkmark

Permission Group	Permission Name	Lab Mgr	Lab Tech	Scientist
Document Editing	Edit Reader Settings			\checkmark
	Edit Notes Text		~	\checkmark
	Read Empty Plates/Cuvettes		~	\checkmark
	Function Editor			\checkmark
	Overwrite Plate/Cuvette Data			\checkmark
	Edit Graphs, Summaries, and Reductions			~
	Assign Plate Layouts		~	\checkmark
	Edit Sample and Group Information		~	~
	Add/Delete Groups			\checkmark
	Edit Formulas			\checkmark
	Edit Print Options			\checkmark
	Lock/Unlock Sections			\checkmark
	Mask/Unmask Wells		~	\checkmark
	Create/Save As Data Document	~		\checkmark
	Create/Save As Protocol			\checkmark
	Change Auto Save Settings			\checkmark
	Change Auto Export Settings			\checkmark
Statements Signatures	Add/Modify Statements			\checkmark
	Sign Statements	~	~	\checkmark
	Revoke Own Signature	~	~	\checkmark
	Revoke Any Signature			
Document Management	Rename Document	\checkmark		\checkmark
	Move Document	~		\checkmark
	Delete Document	~		\checkmark
	Unlock Document	\checkmark		
	Import Documents			\checkmark
	Export Documents	~	~	\checkmark

Default Role Permissions (continued)

Permission Group	Permission Name	Lab Mgr	Lab Tech	Scientist
Folder Management	Add Folder	~		\checkmark
	Rename Folder	~		\checkmark
	Move Folder	~		\checkmark
	Delete Empty Folder	~		\checkmark
	Hide/Unhide Folder	~		\checkmark
Instrument	Lock/Unlock Instrument	~		\checkmark
	Instrument Simulator			\checkmark

Default Role Permissions (continued)

Appendix B: Audit Events

The GxP Admin Software and the SoftMax Pro GxP Software generate audit trails that record the events that user actions trigger. For each event instance, the audit trail records the date and time, the user name of the person who was logged on, and the event action. Users cannot alter the audit trail, nor can they edit or delete audit trail entries.

Note: StakMax specific commands and scripts you use in the StakMax Software are not tracked in the audit trail and do not have user permission settings. Please implement standard work as required by your organization.

The following events trigger a System Audit Trail entry:

- Application Close on Lock: Application closed while in locked state.
- Auto Export Failed: Auto Export to location failed.
- Auto Export Settings Edited: User changed Auto Export settings.
- Auto Export Started: Auto Export started to export to a single location.
- Auto Export Started with Multiple Locations: Auto Export started to export to multiple locations.
- Auto Read Settings Confirmed: User confirmed Auto Read settings.
- Auto Save Settings Edited: User changed Auto Save settings.
- Cuvette Set Section Added: User added a Cuvette Set section to an experiment.
- Cuvette Set Section Appended Data: User appended data to a Cuvette Set section.
- Cuvette Set Section Data Pasted: User pasted data into a Cuvette Set section.
- Cuvette Set Section Data Replaced: User replaced data in a Cuvette Set section.
- Cuvette Set Section Reference Read: User started a Cuvette Set reference read.
- Data Section Data Deleted: User deleted data from a Plate section or a Cuvette Set section.
- Data Section Data Pasted Over Existing: User pasted data over existing data in a Plate section or a Cuvette Set section.
- **Document Unlocked**: User unlocked a document.
- Document Altered, Modified, or Tampered: Document has been changed, modified, or tampered with outside of the SoftMax Pro Software.
- **Document Audit Trail Import Error**: An error occurred importing a document audit trail.
- Document Audit Trail Note Added: User added a not to the audit trail.
- **Document Changes Discarded**: User discarded changes in the document.
- **Document Closed**: User closed a document.
- **Document Created**: User created a new document.
- Document Created from Non-Compliant: Document was created from a noncompliant document.

- **Document Deleted**: User deleted a document from the database.
- **Document Exported**: User exported a document in SoftMax Pro format.
- **Document Legacy Audit Trail Imported**: The audit trail from an imported document was successfully imported to the database.
- **Document Legacy Imported**: Documents were successfully imported into the database.
- **Document Moved**: User moved a document to a different location in the database.
- **Document Open Failed**: An error occurred when a user attempted to open a document.
- Document Opened: User opened a document.
- **Document Print Canceled**: User canceled the print of a document.
- **Document Printed**: User printed a document.
- Document Protocol Created: User created a new protocol.
- ***Document Released Protocol Created**: User created a Released protocol from a Reviewed document.
- **Document Renamed**: User changed the name of a document.
- **Document Save**: User saved a document using Save or Save As.
- **Document Save As PDF**: User saved a document in a .pdf format.
- **Document Save Failed**: An error occurred when a user attempted to save a document.
- Document Selected Sections Exported: User exported selected sections.
- ***Document Status Changed**: User changed the status of a document.
- **Document Unlocked**: User unlocked a document.
- **Document Workflow Activity**: Workflow activity in a document captured.
- Experiment Added: User added an experiment to a document.
- Experiment Deleted: User deleted an experiment from a document.
- **Experiment Duplicated**: User duplicated an experiment in a document.
- Experiment Read Started: User started a read.
- **Experiment Renamed**: User changed the name of an experiment in a document.
- Experiment Section Deleted: User deleted a section from an experiment.
- **Experiment Section Duplicated**: User duplicated a section in an experiment.
- Experiment Section Locked: User locked a section in an experiment.
- Experiment Section Print Disabled: User marked a section as not printable.
- Experiment Section Print Enabled: User marked a section as printable.
- **Experiment Section Renamed**: User changed the name of a section in an experiment.
- Experiment Section Unlocked: User unlocked a section in an experiment.
- **Experiment Summary Formula Added or Edited**: User added or edited a Summary formula in a Group section or a Note section.
- Experiment Summary Formula Deleted: User deleted a Summary formula from a Note section or a Group section.

- Experiment Summary Formula Pasted: User pasted a Summary formula into a section in an experiment.
- Folder Added: User added a folder to the database.
- Folder Deleted: User deleted a folder from the database.
- Folder Hidden: User hid a folder in the database.
- Folder Moved: User moved a folder to a new location in the database.
- Folder Renamed: User changed the name of a folder in the database.
- Folder Unhidden: User unhid a folder in the database.
- **Graph Section Added**: User added a Graph section to an experiment.
- Graph Section Curve Fit Changed: User changed the curve fit in a Graph section.
- Graph Section Settings Edited: User changed the settings in a Graph section.
- **Group Section Column Formula Added or Edited**: User added or edited a Column formula in a Group section.
- Group Section Column Pasted: User pasted a column into a Group section.
- Group Section Cut or Deleted Column: User removed a column from a Group section.
- Group Section Settings Edited: User changed the settings in a Group section.
- Instrument Calibrated: User calibrated an instrument.
- Instrument Locked: User locked the touchscreen on a SpectraMax iD3 or a SpectraMax iD5.
- Instrument Selected: User selected an instrument to acquire data.
- Instrument Temperature Turned Off: User turned the instrument temperature control off.
- Instrument Temperature Turned On: User turned the instrument temperature control on.
- Note Section Added: User added a Note section to an experiment.
- Plate Section Added: User added a Plate section to an experiment.
- Plate Section Append Read: User started a read to append data.
- Plate Section Append Data: User appended data to a Plate section.
- Plate Section Clone Added: User cloned a plate in a Plate section.
- Plate Section Clone Delete Confirmed: User confirmed the deletion of a plate clone.
- Plate Section Clone Deleted: User deleted a plate clone from a Plate section.
- Plate Section Data Import Failed: Data import into a Plate section failed.
- Plate Section Data Import Notification: Data imported into a Plate section contains saturated or invalid values.
- Plate Section Data Imported: Data was imported into a Plate section.
- Plate Section Data Pasted: User pasted data into a Plate section.
- Plate Section Data Replaced: User replaced data in a Plate section.
- Plate Section Read: User started a read.
- Plate Section Read Canceled: User canceled a read.

- Plate Section Read Finished: Read finished.
- Plate Section Read Interrupted: User interrupted a read.
- Plate Section Reduction Changes Accepted: User confirmed changes made to reduction settings.
- Plate Section Settings Edited: User changed the settings for a Plate section.
- Plate Section Template Edited: User changed a plate template.
- Plate Section Template Imported: User imported a template into a Plate section.
- Plate Section Template Pasted: User pasted a template into a Plate section.
- Plate Section Wells Masked: User masked wells in a Plate section.
- Plate Section Wells Unmasked: User unmasked wells in a Plate section.
- Print Options Edited: User changed print options.
- **Project Activated**: User activated a Project.
- **Project Added**: User added a Project.
- **Project Deactivated**: User deactivated a Project.
- ***Project Edited**: User changed a Project.
- Role Activated: User activated a Role.
- Role Added: User added a new Role.
- Role Deactivated: User deactivated a Role.
- *Role Edited: User changed a Role.
- Statement Added: User added a statement to a document.
- Statement All Signatures Removed: All signatures removed when a new document was created by Save As.
- Statement Deleted: User deleted a statement from a document.
- Statement Edited: User edited a statement.
- Statement Pasted: User pasted a statement into a document.
- **Statement Revoke Signature Canceled**: User did not enter their password when revoking a signature from a statement and the signature was not revoked.
- Statement Signed: User signed a statement.
- **Statement Signed Statement Canceled**: User did not enter their password when adding a signature to a statement and the statement was not signed.
- **Statement Signed Statement Removed**: Signed statement removed in the process of creating a new document from a released protocol.
- Statement Single Signature Revoked: User revoked a single signature from a statement.
- System Database Restored: User restored the database.
- **System EDB File Imported**: User imported a file that contains the usernames of users from legacy GxP Admin Software.
- System Offline Data Sync Ended: The SoftMax Pro Software finished synchronizing offline data.
- System Offline Status: The SoftMax Pro software went into offline status.

- ***System Server License Installed**: User added additional SoftMax Pro Software access licenses.
- ***System Settings Changed**: User changed settings in the GxP Admin Portal.
- User Activated: User activated a user.
- User Added: User added a new user.
- ***User Admin Permission Granted**: User was granted the GxP Admin Access permission.
- ***User Admin Permission Revoked**: User revoked the GxP Admin Access permission from a user.
- User Blocked: User blocked due to too many incorrect login attempts.
- User Change Password Attempted: User attempted to change their password.
- User Change Password Canceled: User started to change their password but did not complete the process.
- User Change Password Successful: User successfully changed their password.
- User Deactivated: User deactivated a user.
- ***User Edited**: User changed a user.
- User License Assigned: User was assigned a license to access the SoftMax Pro software.
- User License Unassigned: User removed a SoftMax Pro access license from a user.
- User Logged In: User has logged into the GxP Admin Portal or the SoftMax Pro Software.
- User Logged Out: User has logged out.
- User Logon Attempted: User tried to log into the software.
- ***User Logon Failed**: User unsuccessfully attempted to log into the GxP Admin Portal or the SoftMax Pro Software.
- User Logon Successful: User successfully logged into the software.
- User Password Changed: User changed their password.
- User Software Unlock Failed: An error occurred when a user attempted to unlock the software.
- User Software Unlocked: User unlocked the software.
- * Can trigger a notification for the Home page.

Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

Our Support website, www.moleculardevices.com/service-support, has a link to the Knowledge Base, which contains technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance after consulting the Knowledge Base, you can submit a request to Molecular Devices Technical Support.

You can contact your local representative or Molecular Devices Technical Support at 800-635-5577 X 1815 (North America only) or +1 408-747-1700. In Europe call +44 (0) 118 944 8000.

To find regional support contact information, visit www.moleculardevices.com/contact.

Contact Us

Phone: +1-800-635-5577 Web: moleculardevices.com Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners. Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES ©2020 Molecular Devices, LLC. All rights reserved. 0112-0123 J

