

SoftMax Pro 7.3 GxP – Functional Requirements Specification Assessment

This Functional Requirements Specification document is designed retrospectively to address SoftMax® Pro 7.3 GxP Data Acquisition and Analysis Software as suitable compliance software.

Molecular Devices is an ISO registered company certified to ISO 9001:2015 under registrar BSI certificate number FS 534246.

We provide SoftMax Pro GxP Software that extends Molecular Devices' leading data acquisition and analysis solution into regulated laboratories working under GMP, GLP, 21 CFR Part 11, and other similar guidelines for secure electronic records.

SoftMax Pro GxP Software provides the functionality that enables compliance. SoftMax Pro GxP Software is only a tool which assists customers in becoming 21 CFR Part 11 compliant. It is the responsibility of the customer to comply with 21 CFR Part 11.

Tim Bolus, MPM, MBA
Compliance Product & Program Manager

Shreen Ibrahim
Director, Global Quality

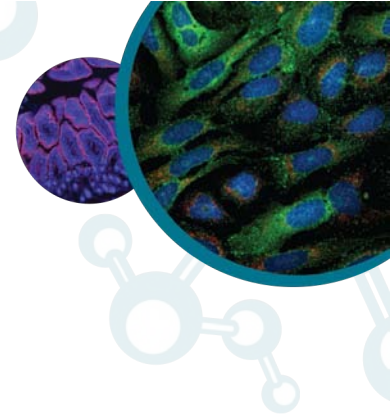
Revision History

1	Original document
2	Added the General category.
3	<ul style="list-style-type: none"> Modified the Curve Fitting section and renamed it as Interpolation & Curve Fitting Removed the Statistical Analysis of Plate Data and Statistical Analysis of Cuvette Set Data sections. Added new items: GEN-03 and GEN-04 Added the Optical Configuration Detection section.

3860 N First Street
San Jose, CA 95134
United States

toll-free 1.800.635.5577
office 408.747.1700

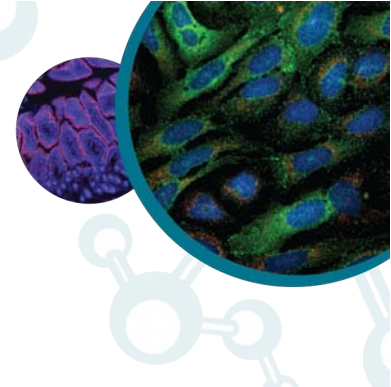
moleculardevices.com



GxP ADMIN PORTAL FUNCTIONALITY		
ID	Requirement	Comments
GXP-01	The software shall permit adjustment of default password settings in GxP Admin Portal.	Permissions can be configured to end-user business processes in GxP Admin Portal.
GXP-02	The software shall permit functionality to add users.	
GXP-03	The software shall permit functionality to edit users.	
GXP-04	The software shall permit access to the GxP Admin Portal.	
GXP-05	The software shall permit changing the GxP Admin Portal first login password.	
GXP-06	The software shall permit modifying the date and time format.	
GXP-07	The software shall permit access to the System Audit Trail.	
GXP-08	The software shall permit assigning access to the GxP Admin Portal.	
GXP-09	The software shall permit changing a user password in the GxP Admin Portal.	
GXP-10	The software shall permit functionality to add Roles.	
GXP-11	The software shall permit functionality to edit Roles.	
GXP-12	The software shall permit functionality to copy Roles.	
GXP-13	The software shall permit functionality to add Projects.	
GXP-14	The software shall permit functionality to edit Projects.	
GXP-15	The software shall permit functionality to copy Projects.	
GXP-16	The software shall provide shortcuts to the Home page.	
GXP-17	The software shall permit access to the SoftMax Pro GxP software.	
GXP-18	The software shall provide Administrator Contact information.	
GXP-19	The software shall permit functionality to automatically lock out users after a defined period of inactivity.	
GXP-20	The software shall permit a maximum number of false login attempts.	
GXP-21	The software shall allow management of notifications.	

CUSTOM CREDENTIALS		
ID	Requirement	Comments
CCP-01	The software shall allow a minimal length of characters for passwords.	Permissions can be configured to end-user business processes in GxP Admin Portal.
CCP-02	The software shall allow uppercase letters for passwords.	
CCP-03	The software shall allow lowercase letters for passwords.	
CCP-04	The software shall allow digits for passwords.	
CCP-05	The software shall allow special characters for passwords.	
CCP-06	The software shall allow that passwords differ from username.	
CCP-07	The software shall maintain a history of previous passwords.	
CCP-08	The software shall allow password aging.	

WINDOWS ACTIVE DIRECTORY CREDENTIALS		
ID	Requirement	Comments
WC-01	The software shall allow use of Windows Active Directory credentialing.	Permissions can be configured to end-user business processes in GxP Admin Portal.
WC-02	The software shall allow use of Windows Active Directory password.	

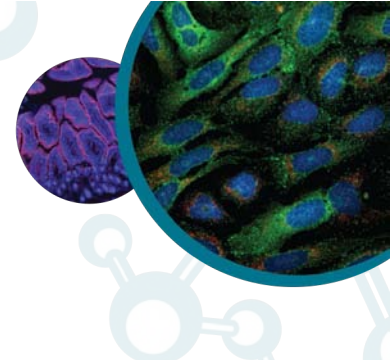


FOLDER MANAGEMENT		
ID	Requirement	Comments
FM-01	The software shall allow for creation of folders.	Permissions can be configured to end-user business processes in GxP Admin Portal.
FM-02	The software shall allow for movement of folders.	
FM-03	The software shall allow for the hiding/unhiding of folders.	
FM-04	The software shall allow for renaming of folders.	
FM-05	The software shall allow for deletion of empty folders.	

BUILD DOCUMENT WORKFLOW		
ID	Requirement	Comments
BDW-01	The software shall allow users to create/save data document(s).	Permissions can be configured to end-user business processes in GxP Admin Portal.
BDW-02	The software shall allow users to create/save protocol document(s).	
BDW-03	The software shall have auto-saving capability.	
BDW-04	The software shall have auto-export capability.	

DOCUMENT MANAGEMENT WORKFLOW		
ID	Requirement	Comments
DMW-01	The software shall allow movement of documents within the project folders.	Permissions can be configured to end-user business processes in GxP Admin Portal.
DMW-02	The software shall allow for renaming of documents within the project folders.	
DMW-03	The software shall allow the deletion of documents within the project folders.	
DMW-04	The software shall allow the import of documents into project folders.	
DMW-05	The software shall allow the export of documents from project folders.	
DMW-06	The software shall allow the ability to open imported documents in the project folders.	

DOCUMENT EDITING WORKFLOW		
ID	Requirement	Comments
DEW-01	The software shall allow a user to open a document that has the status: 'In Work'	Permissions can be configured to end-user business processes in GxP Admin Portal.
DEW-02	The software shall allow a user to edit plate reader settings.	
DEW-03	The software shall allow a user to edit graphs, summaries and reductions.	
DEW-04	The software shall allow a user to edit Note section text	
DEW-05	The software shall allow a user to simulate raw data acquisition.	
DEW-06	The software shall allow a user to read empty plates or cuvette sets.	
DEW-07	The software shall allow a user to assign plate layouts.	
DEW-08	The software shall allow a user to edit Sample and Group information.	
DEW-09	The software shall allow a user to add/delete assigned Groups.	
DEW-10	The software shall allow a user to edit formulas.	
DEW-11	The software shall allow a user to lock/unlock sections of a document.	
DEW-12	The software shall allow a user to add/modify e-signature statements.	
DEW-13	The software shall allow a user to edit printing options.	
DEW-14	The software shall allow a user to overwrite plate or cuvette set data.	
DEW-15	The software shall allow a user to mask/unmask wells in a Plate section of a document.	



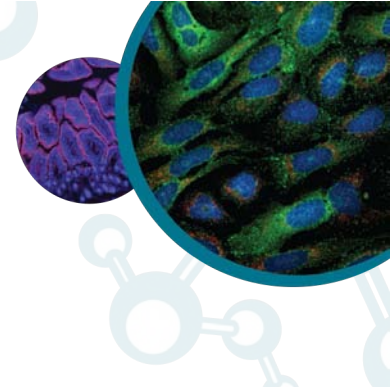
STATEMENTS & SIGNATURES WORKFLOW		
ID	Requirement	Comments
SSW-01	The software shall allow users to sign e-signature statements.	Permissions can be configured to end-user business processes in GxP Admin Portal.
SSW-02	The software shall allow users to revoke any e-signature statement.	
SSW-03	The software shall allow users to revoke their own e-signature.	

DOCUMENT REVIEW WORKFLOW		
ID	Requirement	Comments
DRW-01	The software shall allow a user to set a document to the status: 'Review Pending'.	Permissions can be configured to end-user business processes in GxP Admin Portal.
DRW-02	The software shall allow a user to open a document that has the status: 'Review Pending'.	
DRW-03	The software shall allow a user to unlock a document.	
DRW-04	The software shall allow a user to set a document to the status: 'Reviewed'.	
DRW-05	The software shall allow a user to open a document that has the status: 'Reviewed'.	
DRW-06	The software shall allow a user to set a document to the status: 'Canceled'.	
DRW-07	The software shall allow a user to open a document that has the status: 'Canceled'.	
DRW-08	The software shall allow a user to release a protocol document.	

GENERATE DATA WORKFLOW		
ID	Requirement	Comments
GDW-01	The software shall allow a user to open a document that has the status: 'Released'.	Permissions can be configured to end-user business processes in GxP Admin Portal.
GDW-02	The software shall allow a user to open a document that has the status: 'Open In Progress'.	
GDW-03	The software shall allow a user to generate compliance data.	

DOCUMENT APPROVAL WORKFLOW		
ID	Requirement	Comments
DAW-01	The software shall allow a user to set a document to the status: 'Set Approval Pending'.	Permissions can be configured to end-user business processes in GxP Admin Portal.
DAW-02	The software shall allow a user to open a document that has the status: 'Set Approval Pending'.	
DAW-03	The software shall allow a user to set a document to the status: 'Approved'.	
DAW-04	The software shall allow a user to open a document that has the status: 'Approved'.	
DAW-05	The software shall allow a user to set a document to the status: 'Outdated'.	
DAW-06	The software shall allow a user to open a document that has the status: 'Outdated'.	
DAW-07	The software shall allow a user to send a document back to the status: 'In Progress'.	

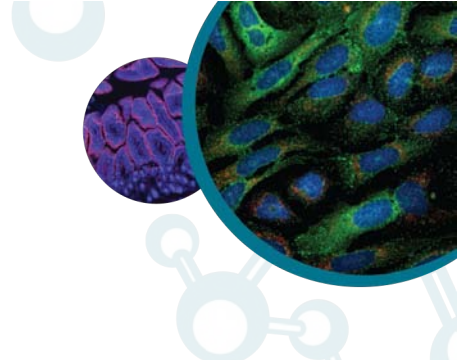
GENERAL		
ID	Requirement	Comments
GEN-01	The software shall suspend calculations and recalculate values.	Permissions can be configured to end-user business processes in GxP Admin Portal.
GEN-02	The software shall allow a user to create custom calculations.	
GEN-03	The software shall collect data in endpoint, kinetic, spectrum, and well scan modes in the automated workflow feature.	
GEN-04	The software shall permit the export of bulk data files outside the database.	



CONNECTIVITY	
<i>Connectivity between the SoftMax Pro GxP Software Suite is out of the box functionality.</i>	
ID	Requirement
CON-01	The SoftMax Pro GxP software can still acquire data when GxP Admin (services) is offline.
CON-02	There shall be connectivity between SoftMax Pro GxP software and a microplate reader.
CON-03	The software shall provide automatic data recovery.
CON-04	The software shall implement the Unlock/Lock instrument permissions to disable the touchscreen on the SpectraMax iD3 or SpectraMax iD5.

INTERPOLATION & CURVE FITTING		
<i>Interpolation and Curve fitting are out of the box functionality of the SoftMax Pro software.</i>		
ID	Requirement	Equation
CFT-01	The software shall interpolate values for a Linear curve fit.	Where $y = A + Bx$
CFT-02	The software shall interpolate values for a Cubic curve fit.	Where $y = A + Bx + Cx^2 + Dx^3$
CFT-03	The software shall interpolate values for a Quadratic curve fit.	Where $y = A + Bx + Cx^2$
CFT-04	The software shall interpolate values for a Quartic curve fit.	Where $y = A + Bx + Cx^2 + Dx^3 + Gx^4$
CFT-05	The software shall interpolate values for a Log-Log curve fit.	Where $\text{Log}(y) = A + B \cdot \text{Log}(X)$
CFT-06	The software shall interpolate values for a Semi-Log curve fit.	Where $y = A + B \cdot \text{Log}(X)$
CFT-07	The software shall interpolate values for a Power curve fit.	Where $y = A + Bx^C$
CFT-08	The software shall interpolate values for a Michaelis-Menten curve fit.	Where $y = Ax/x+B$
CFT-09	The software shall allow for a Log-Logit curve fit.	Where $y = D + [(A-D) / (1 + (x/C)^B)]$
CFT-10	The software shall interpolate values for a 4-Parameter curve fit.	
CFT-11	The software shall interpolate values for a 4-Parameter Global Fit (PLA) curve fit.	
CFT-12	The software shall allow for a 4-Parameter Global Fit (Shared) curve fit.	
CFT-13	The software shall allow for a 5-Parameter curve fit.	Where $y = D + [(A-D) / \{1 + (x/C)^B\}^G]$
CFT-14	The software shall allow for a 5-Parameter Global Fit (PLA) curve fit.	
CFT-15	The software shall allow for a 5-Parameter Global Fit (Shared) curve.	

OPTICAL CONFIGURATION DETECTION	
<i>Optical Configuration for detection cartridges is out of the box functionality of the SoftMax Pro software.</i>	
ID	Requirement
OCD-01	The software shall allow the optical configuration for the Multi-Mode detection cartridge(s).
OCD-02	The software shall allow the optical configuration for the AlphaScreen detection cartridge(s).
OCD-03	The software shall allow the optical configuration for the Fluorescence Intensity detection cartridge(s).
OCD-04	The software shall allow the optical configuration for the Fluorescence Polarization detection cartridge(s).
OCD-05	The software shall allow the optical configuration for the Glow Luminescence detection cartridge(s).
OCD-06	The software shall allow the optical configuration for the Time-Resolved Fluorescence detection cartridge(s).
OCD-07	The software shall allow the optical configuration for the ScanLater detection cartridge(s).

**GXP ADMIN BACKUP**

Settings can be configured to end-user business processes in the GxP Admin Backup tool. Applicable only for stand-alone (single PC) systems.

ID	Requirement
DBB-01	There shall be functionality to backup the GxP Admin Software database.
DBB-02	There shall be functionality to schedule a GxP Admin Software database backup.
DBB-03	There shall be functionality to restore a GxP Admin database.