



SoftMax Pro GxP

SoftMax Pro 7.1 GxP Data Acquisition and Analysis Software

Meet FDA guidelines in GMP/GLP labs with complete validation tools

KEY BENEFITS

- Track and record all changes with the system audit trail
- Enable enterprise-level document sharing with Microsoft SQL database
- Gain complete control over file access permissions
- Maintain data integrity with eSignatures and document workflows
- Reduce IT support with Windows active directory

SoftMax Pro 7.1 GxP Software

SoftMax® Pro 7.1 GxP Software is the latest, most secure software to achieve full FDA 21 CFR Part 11 compliance with streamlined workflows to ensure data integrity. Every step is optimized to simplify analysis and reporting to support our microplate readers.

Our expert team will partner with you to set up single- or enterprise-level software, and provide IQOQ services using our validation package to establish full compliance on your microplate readers. Major data privacy and security improvements support the latest GDPR regulations.

GxP tools for compliance

FDA 21 CFR Part 11 compliance tools are provided for both data acquisition and analysis for 20+ Molecular Devices microplate readers.



Compliance tools

- System audit trail
- Secure data storage in a Microsoft SQL database
- Granular level of control over user permissions
- eSignatures and time stamps



IQ/OQ/PM

- Installation Qualification (IQ) verifies proper installation and connection to SoftMax Pro GxP Software
- Operational Qualification (OQ) tests the mechanical, electrical, and optical functions
- Preventive Maintenance (PM) ensures best working conditions for each instrument



Software validation

SoftMax Pro GxP Validation Package provides comprehensive documentation and tools to validate the software and GxP Admin features.



- Validation time is shortened from 6 months to 3 days using an electronic manual
- Comprehensive tools confirm calculations, curve fits, and parallel line analysis
- Sample data sets shorten authentication and analysis time

SpectraTest Validation Plates

The SpectraTest® Validation Plates provide automated, comprehensive, and traceable validation of microplate reader performance for absorbance, fluorescence, and luminescence read modes.

[The Multi-Mode Validation Plate offers additional testing of TRF, HTRF, and fluorescence polarization read modes on FilterMax™ F3/F5 and SpectraMax® i3x, iD5 and Paradigm® microplate readers.]



Automated for ease of use

All test measurements and calculations are handled automatically by the SoftMax Pro GxP Software protocols



NIST traceability

Calibration of a plate's filter standards is accomplished through the use of an instrument calibrated with primary NIST standards



Recertification service

Validation plates are available for recertification according to ISO 17025 guidelines

Data acquisition and analysis



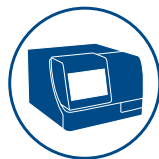
- View data in numeric or scientific notation format, 3D bar graph, heat map, or gray scale modes
- Analyze data using advanced curve-fitting capabilities, relative potency and IC_{50} determination, Michaelis-Menten enzyme kinetics, area under the curve calculations, V_{max} , and more
- Test for outliers using Rosner (ESD) and ROUT methods
- Validate hardware quickly and easily with built-in validation protocols for primary detection modes

Professional services



- MDCare GxP Support Plan for advanced troubleshooting analysis
- Virtual installation services
- Customized protocol support

Support for all our microplate readers



- SpectraMax® iD3/iD5, i3x, M Series, Multi-Mode Microplate Readers
- SpectraMax® L Microplate Reader
- FlexStation® 3 Multi-Mode Microplate Reader
- SpectraMax® ABS and ABS Plus Microplate Readers
- And more

Product	Non-expiring license
SoftMax Pro GxP Software Windows 7/8/10 compatible	<p>Latest version of SoftMax Pro 7 GxP Software, 3 software installations for every purchased user license, GxP Admin Software suite, validation pack, user license certificate, compliance certificate</p> <p>Software part number: SMP7X GXP*</p> <p>User license part number: SMP7 GXP License (3-10) SMP7 GXP License (11-50) SMP7 GXP License (>50)</p> <p>*Requires purchase of any amount of user licenses (minimum three). Tiered pricing is based on number of user licenses purchased. Please select one user license part number when ordering.</p>

Validation	Part number
SpectraTest ABS1 Absorbance Validation Plate	0200-6117
SpectraTest FL1 Fluorescence Validation Plate	0200-5060
SpectraTest LM1 Luminescence Validation Plate	0200-6186
Multi-Mode Validation Plate	0200-7200
Cuvette Validation Set	0200-2420

Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1.800.635.5577	China (Beijing) +86.10.6410.8669	Japan (Osaka) +81.6.7174.8331
United Kingdom +44.118.944.8000	China (Shanghai) +86.21.3372.1088	Japan (Tokyo) +81.3.6362.5260
Europe* 00800.665.32860	Hong Kong +852.3971.3530	South Korea +82.2.3471.9531

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland