



SoftMax Pro 7 GxP Software

KEY FEATURES

- **Validation package minimizes the time and cost of software validation**
- **Admin module controls end user access and usage**
- **Built-in features allow access by auditors without the need to purchase additional licenses**

Trusted

- SoftMax® Pro 7 GxP is our fourth generation software with compliance tools
- Over 100 field and internal staff with full understanding of hardware, applications, and analysis with over a century of combined experience
- More than 160,000 software licenses sold since 2004
- Satisfied customers include all 50 of the top 50 global pharmaceutical companies
- Support from a single vendor that develops both hardware and software offering over 160 specialized protocols for analyzing microplate data

Save time and reduce cost

- Extensive suite of tools can reduce the cost and time of validation by 50% as compared to using multiple platforms to collect and analyze data
- Combines end-to-end chain of custody from capture through analysis to validation of data all on one platform
- Eliminates extra time and loss of accuracy associated with transfer between software platforms
- Robust turnkey solution saves up to \$200K in method development cost vs. alternatives

Updated compliance-focused software

Time and money saving GxP features	2003 (v4)	2005 (v5)	2012 (v6)	2016 (v7)
Audit Trail	✓	✓✓	✓✓	✓✓
Password aging	✓	✓	✓	✓
Granular permission levels	✓	✓✓	✓✓✓	✓✓✓
Curve fitting	6	9	21*	21*
Parallel line analysis (PLA) and relative potency		Protocols	Built-in validation	Built-in validation
Designate secondary administrators		✓	✓	✓
Software validation package	✓	✓	✓✓	✓✓
Number of readers supported	14	18	22	23
FlexStation 3 support		✓		✓
Operating systems supported	Win XP	Win XP/7	Win 7/8	Win 7/8/10
IQ/OQ	Paper	Paper	Digital	Digital
Confidence interval calculations			✓	✓
Parameter dependency to determine best curve fit			✓	✓
Enhanced note section with text placement flexibility			✓	✓
Customized reporting with images, live graphs, and tables			✓	✓
Print preview and custom header and footer options			✓	✓✓
Context-sensitive searchable help			✓	✓
Import file format for data analysis	.txt	.txt	.txt, .xml	.txt, .xml
Export file format for transport to database	.txt	.txt, .xml	.txt, .xls, .xml	.txt, .xls, .xml
Multitask kinetics with execution log				✓
Validation plate protocols for i3(x) and Paradigm cartridges and FilterMax F3/F5				✓

*Enhanced graphical options and 21 curve fits including 4- and 5-parameter logistic models

Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Data acquisition and analysis compliance software

- 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
- Validate hardware quickly and easily with built-in validation protocols for primary detection modes
- Sign password-protected electronic signatures for data file verification, authorization, and approval
- Document the history of user actions for each data file in computer-generated, time-stamped audit trails
- Allow remote user access to SoftMax Pro 7 GxP Software through the TCP/IP connection option
- Take advantage of our expert installation and custom protocol services to speed up deployment and adoption

Professional Services

MDCare GxP Support Plan

- Priority response
- Advanced analysis support
- Up to 2 days of lab support (Molecular Devices site only)
- Remote instrument access
- User accounts and permissions troubleshooting
- Advanced support specifically for compliant labs with validation questions on SoftMax Pro and instruments

Virtual Installation Services

- Feedback on customer IT system architectures including virtualization
- Installation issue troubleshooting
- MDC file server licensing setup and configuration
- GxP Admin permission and user account setup help
- Licensing setup and configuration

Custom Protocol Services

- Convert current protocols to SoftMax Pro or develop new protocols
- Use our Ph.D. level experts to create SoftMax Pro protocols suited for your exact needs
- Standard, mid-level, and high-level complexity protocol creation options