

## DATA SHEET

# SoftMax Pro 7.1.1 GxP Software

Acquisition and analysis software for microplate readers  
in compliance labs



SoftMax  
Pro

## Trusted

- SoftMax® Pro 7.1.1 GxP is our latest generation software with compliance tools
- Satisfied customers include all 50 of the top 50 global pharmaceutical companies
- Our global field and development teams have broad expertise in laboratory equipment, applications, and analysis
- More than 160,000 software licenses sold since 2004
- 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

## KEY FEATURES

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

## Updated compliance-focused software

Time and money saving GxP features	2005 (v5.4)	2012 (v6.5.1)	2016 (7.0.3)	2020 (v7.1.1)
System audit trail (Event tracking with date and time stamps)				✓
Secure file database (SQL)				✓
Administrative control of granular permission levels	User based	User based	User based	Role based
Integrated password settings (LDAP)				✓
Curve fitting options	9	21*	21*	22**
Integrated tools for parallel line analysis (PLA) and relative potency	✓	✓✓	✓✓	✓✓✓
Import file format for data analysis	.txt	.txt, .xml	.txt, .xml	.txt, .xml
Export file format for transport to database	.txt, .xml	.txt, .xls, .xml	.txt, .xls, .xml, PDF	.txt, .xls, .xml, PDF
Legacy file support		v5	v5, v6	v6, v7
Operating systems supported	Win XP/7	Win 7/8	Win 7/8/10	Win 10
Server operating systems supported				Win 2016 & 2019
Virtual environment support				✓
Validation plate protocols for i3(x) and Paradigm cartridges and FilterMax F3/F5			✓	✓
SpectraMax® ABS and ABS Plus reader support				✓
SpectraMax® L reader support	✓		✓	✓

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

\*\*Custom curve fit option

## Professional Services

### MDCare GxP Support Plan

- Advanced analysis support
- Remote instrument troubleshooting
- Advanced support specifically for compliant labs with validation questions on SoftMax Pro and instruments
- Priority response

### Virtual Installation Services

- Feedback on customer IT system architectures including virtualization
- Consultation on acceptable IT infrastructure for deployment
- Remote installation support
- Troubleshooting installation issues

### 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

### Contact Us

Phone: +1.800.635.5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto: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