

DATA SHEET

SoftMax Pro 7.1.2 GxP Software

Acquisition and analysis software for microplate readers
in compliance labs



SoftMax
Pro

Trusted

- SoftMax® Pro 7.1.2 GxP Software 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)	2021 (v7.1.2)
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	Role based
Integrated password settings (LDAP)				✓	✓
Curve fitting options	9	21*	21*	22**	22**
Integrated tools for parallel line analysis (PLA) and relative potency	✓	✓✓	✓✓	✓✓✓	✓✓✓
Import file format for data analysis	.txt	.txt, .xml	.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	.txt, .xls, .xml, PDF
Legacy file support ¹		v5	v5, v6	v6, v7	v4, v5, v6, v7
Operating systems supported	Win XP/7	Win 7/8/10	Win 7/8/10	Win 10	Win 10
Server operating systems supported				Win 2016 & 2019	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	✓		✓	✓	✓
Instrument Diagnostics Mode (for SpectraMax® i3x, SpectraMax® iD3, SpectraMax® iD5, SpectraMax® Paradigm)					✓

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

**Custom curve fit option

¹Legacy file formats supported in 7.1.2: SMP 4.0 – 5.4 Data (*.pda); SMP 4.0 – 5.4 Protocol (*.ppr); SMP 4.3 – 5.4 GxP Data (*.edaj); SMP GxP 4.3 – 5.4 Protocol (*.epr)

Professional Services

SMPCare: GxP Support Plan

- Advanced analysis support
- Remote instrument troubleshooting
- Advanced support specifically for compliant labs with validation questions on SoftMax Pro Software
- 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 Software or develop new protocols
- Use our Ph.D. level experts to create SoftMax Pro Software 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
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1.800.635.5577 Taiwan/Hong Kong +886.2.2656.7585
 United Kingdom +44.118.944.8000 Japan +81.3.6362.9109
 Europe* 00800.665.32860 South Korea +82.2.3471.9531
 China +86.4008203586 India +91.73.8661.1198

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