

DATA SHEET

SoftMax Pro GxP Software

Acquisition and analysis software for microplate readers
in compliance labs



SoftMax
Pro

Trusted

- SoftMax® Pro GxP Software is our latest generation software that supports 21 CFR Part 11 and Annex 11 compliance for GMP/GLP regulated labs
- 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 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 | 2012 (v6.5.1) | 2016 (7.0.3) | 2020 (v7.1.1) | 2021 (v7.1.2) | 2023 (v7.2) |
|---|------------------|-----------------------|--------------------------------|--------------------------------|--------------------------------|
| System audit trail (Event tracking with date and time stamps) | | | ✓ | ✓ | ✓ |
| Secure file database (SQL) | | | ✓ | ✓ | ✓ |
| Administrative control of granular permission levels | User based | User based | Role based (52 permissions) | Role based (52 permissions) | Role based (53 permissions) |
| Integrated password settings (LDAP) | | | ✓ | ✓ | ✓ |
| Curve fitting options | 21* | 21* | 22 ^a | 22 ^a | 24 ^{a,b} |
| Integrated tools for parallel line analysis (PLA) and relative potency | ✓✓ | ✓✓ | ✓✓✓ | ✓✓✓ | ✓✓✓ |
| Import file format for data analysis | .txt, .xml | .txt, .xml | .txt, .xml | .txt, .xml | .txt, .xml |
| Export file format for transport to database | .txt, .xls, .xml | .txt, .xls, .xml, PDF | .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 | v4, v5, v6, v7 |
| Operating systems supported | Win 7/8/10 | Win 7/8/10 | Win 10 | Win 10 | Win 10/11 |
| Server operating systems supported | | | Win 2016 & 2019 | 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 | | ✓ | ✓ | ✓ | ✓ |
| SpectraMax [®] Mini reader support | | | | | ✓ |
| Instrument Diagnostics Mode (for SpectraMax [®] i3x, SpectraMax [®] iD3, SpectraMax [®] iD5, SpectraMax [®] Paradigm) | | | | ✓ | ✓ |

^aEnhanced graphical options including 4- and 5-parameter logistic models

^bCustom curve fit option

¹Global curve fit option

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

- Use our Ph.D. level experts to convert protocols to SoftMax Pro documents or create new protocols suited for your business 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, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom