

DATA SHEET

SoftMax Pro Software

Acquisition and analysis software for microplate readers



SoftMax
Pro

Trusted

- SoftMax® Pro Software 7 is our latest generation software and has been continuously improved since launch of version 1 in 1995
- SoftMax Pro Software is the most published microplate reader acquisition and analysis software
- Satisfied customers include ALL of the top 50 global pharmaceutical companies
- Our global field and development teams have broad expertise in laboratory equipment, applications, and analysis

KEY FEATURES

- Over 160 built-in protocols to simplify data acquisition and analysis
- Updated protocols available through customer care portal, [Spectranet](#)
- Advanced reader control: Custom assay workflows, multitask kinetics, and pause and resume kinetic reads with discontinuous kinetics feature
- Supports traditional and unconventional microplate reader assays, including imaging cytometry and western blots
- Flexible calculations with data analysis features such as 24 different curve fit options, simplified cross-plate analysis and interpolation, and automatic calculations of relative potency, EC_{50} , and Z factors
- Spectral optimization wizard to maximize assay signal window (SpectraMax® i3x, iD3, iD5, and Paradigm Multi-Mode Microplate Readers)

Updated microplate reader software

Time and money saving SoftMax Pro features	2012 (v6.5.1)	2016 (7.0.3)	2020 (v7.1.1)	2021 (v7.1.2)	2023 (v7.2)	2024 (v7.3)
Curve fitting options	21*	21*	22 ^a	22 ^a	24 ^{a,b}	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	*.txt, *.xml
Export file format	*.txt, *.xls, *.xml	*.txt, *.xls, *.xml, *PDF	*.txt, *.xls, *.xml, *PDF	*.txt, *.xls, *.xml, *PDF	*.xslt, *.txt, *.xls, *.xml, *PDF	*.xslt, *.txt, *.xls, *.xml, *PDF
Legacy file support ¹	v5	v5, v6	v6, v7	v4, v5, 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	Win 10/11
Graph functionality and display for publications	✓	✓	✓✓	✓✓	✓✓✓***	✓✓✓***
Instrument Diagnostics Mode (for SpectraMax [®] i3x, SpectraMax [®] iD3, SpectraMax [®] iD5, SpectraMax [®] Paradigm)				✓	✓	✓
Bulk export for multiple documents						✓
Tool tips translated to Simplified Chinese						✓

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

^aCustom curve fit option

^bGlobal 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)

***graph annotations

Instruments compatible with the SoftMax Pro 7: SpectraMax[®] ABS, ABS Plus, i3, i3x, iD3, iD5, Paradigm, Mini, VersaMax, Plus384, L, M5/e, M4, M3, M2/e, 340PC 384, 190/EXT, Gemini XPS/EM, FlexStation 3, F5, F3, DTX880, DTX800, Vmax, Emax and StakMax

Instruments not supported by SoftMax Pro 7: SpectraMax[®] 190 Absorbance Microplate Reader (Serial Numbers beginning with Nxxxxx are not supported) and VersaMax Microplate Reader (Serial Numbers beginning with Bxxxxx are not supported)

Professional Services

SoftMax Pro Software

- Integration with leading robotics and LIMS partners enables both data analysis and instrument control in automated environments
- Ph.D. level technical support
- SoftMax Pro specific training courses available upon request

Virtual Installation Services for the GxP edition

- 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

