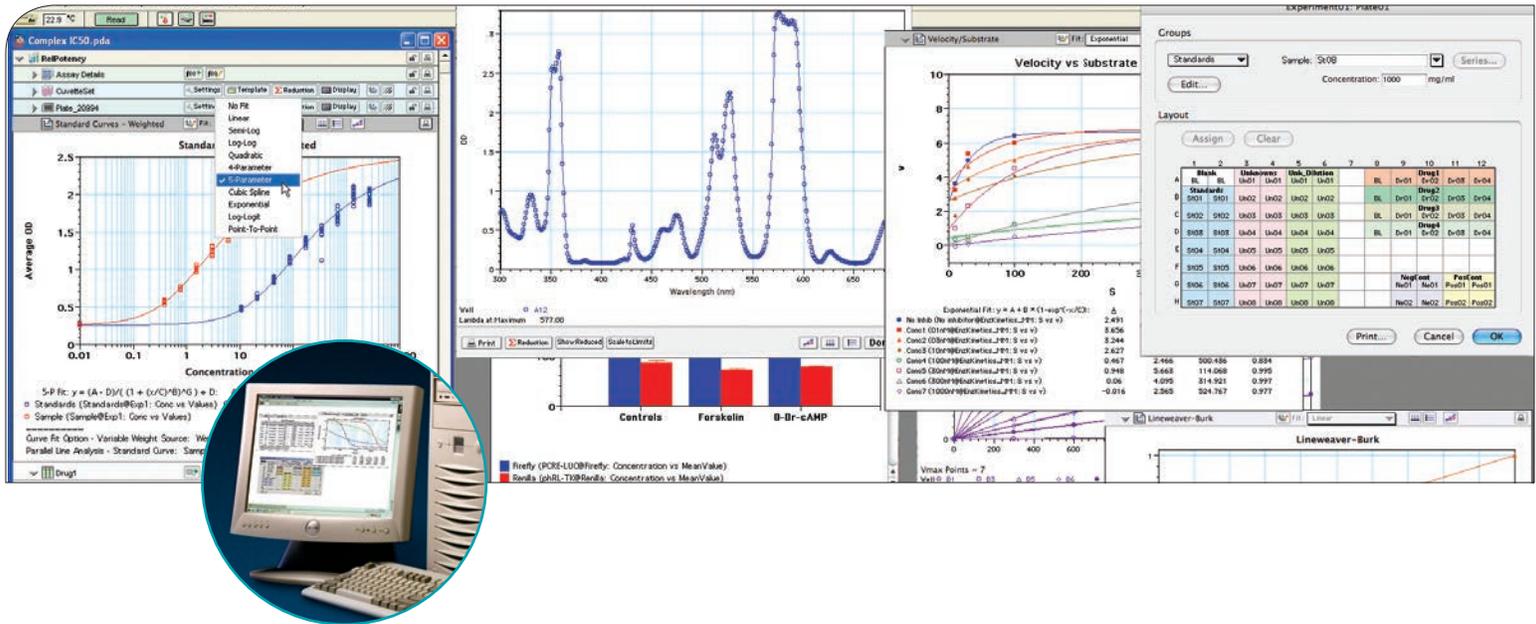


SoftMax Pro 5 Software

THE INDUSTRY STANDARD IN MICROPLATE DATA ANALYSIS



- SIMPLE AND CUSTOMIZABLE ASSAY SETUP
- FLEXIBLE AND POWERFUL DATA ANALYSIS
- COMPLETE FDA 21 CFR PART 11 COMPLIANCE TOOLS
- INTEGRATION CAPABILITIES WITH LEADING ROBOTIC SYSTEMS
- OPTIONAL VALIDATION TOOLS

SoftMax[®] Pro 5 Software from Molecular Devices offers extensive data capture, analysis, and reduction capabilities for its absorbance, luminescence, fluorescence, and multi-detection microplate readers for customers working in pharmaceutical, biotechnology, academia, hospital, or government laboratories.

With over 120 ready-to-run assay protocols for life science research and drug discovery assay development and screening, researchers can customize experiment protocols, data display, analysis, and the formation of meaningful reports. This industry-leading software package can be integrated into Molecular Devices' many partner robotics systems.

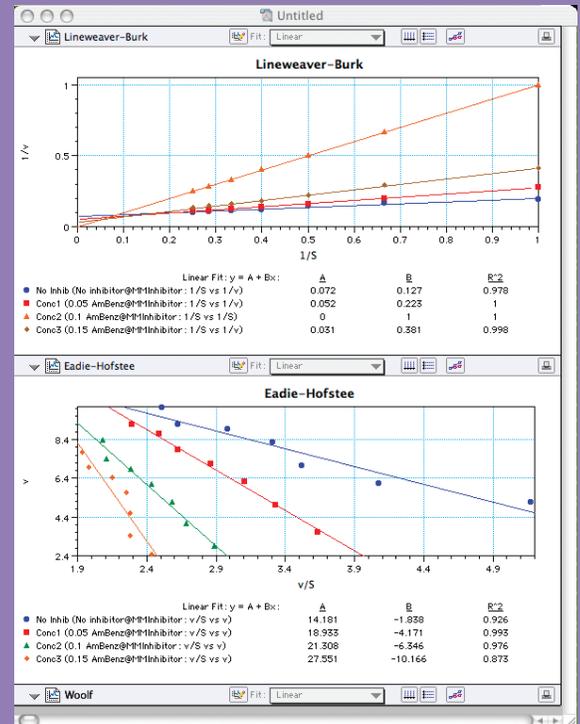
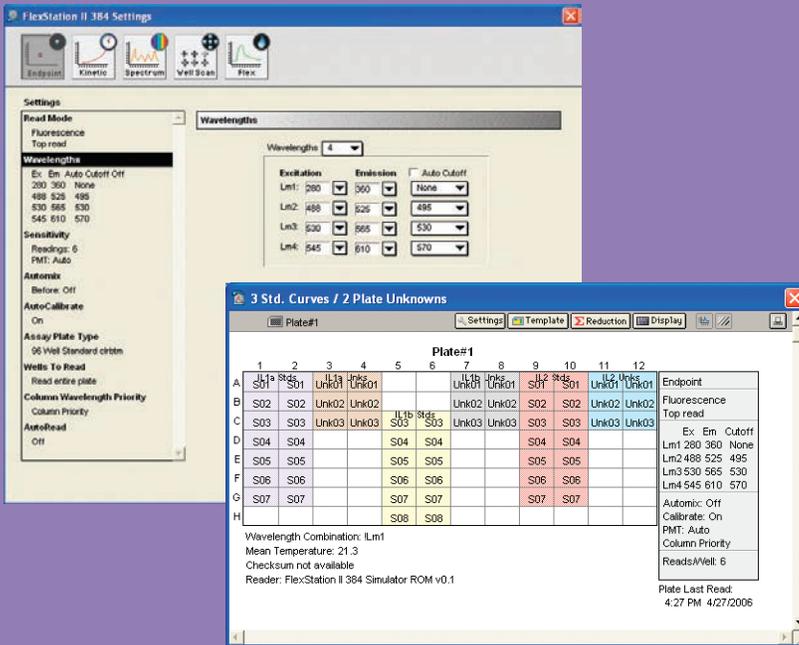
SoftMax Pro 5 Software is truly cross-platform, running on Windows and native to Mac OS X, allowing users to open files interchangeably across platforms.

TWO EDITIONS TO MEET NEEDS

SoftMax Pro 5 Software provides extensive data acquisition and analysis features for Gemini, LMax, SpectraMax[®], VersaMax[™], VMax[®], EMax[®], UVMax, FlexStation[®], and ThermoMax benchtop Microplate Readers. Data collected by Analyst[®] and FLIPR[®] Systems can also be imported and analyzed in the software. SoftMax Pro 5 GxP Software provides all the data collection and analysis tools of the standard edition, along with features critical to operating under GLP/GMP compliance guidelines.

Additional validation tools are available to reduce the time, effort, and cost of certifying Molecular Devices' microplate systems.

Simple and Customizable



Using SoftMax Pro 5 Software is as easy as 1, 2, 3!

1. Turn on the instrument and launch SoftMax Pro 5 Software. The software identifies the instrument and displays only appropriate control options.
2. Open one of the 120+ included protocols and enter template data.
3. Press the Read button, save the file, and print out a report.

The graphing features in SoftMax Pro 5 Software save time and hassle by generating powerful, customized graphs within a single application; create and lay out microplate templates with custom names, descriptions, concentrations, and dilutions.

SIMPLE SETUP

SoftMax Pro 5 Software helps take advantage of an instrument's capabilities by providing:

- Automatic instrument detection
- A wide variety of scan types: Endpoint, Kinetic, Spectral Scanning, and Well Scanning
- The ability to read multiple wavelengths[‡]
- Temperature regulation and patented* AutoMix features to optimize assay environments
- Extensive predefined support for 6- to 384-well plate formats[‡]
- The ability to select specific rows of interest to shorten plate read times
- Patented* PathCheck® technology to automate well volume normalization for OD measurements[§]

CUSTOM TEMPLATES

Create and define plate templates—then view, import, or export them with easy-to-use editing features designed to streamline assay and experimental setup.

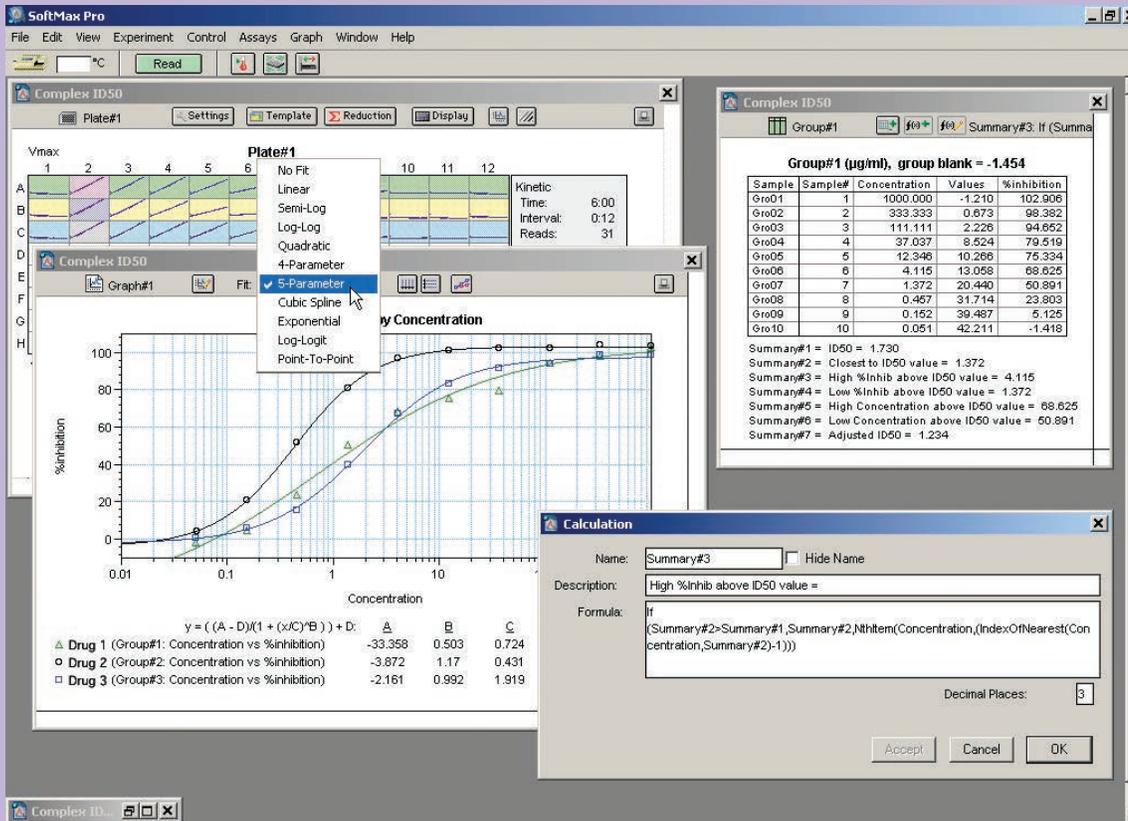
- Assign unilateral samples, standards, controls, or blanks
- Visualize groups quickly with color-coded graphics
- Quickly replicate experiment templates using the Copy/Paste features
- Easily assign dilutions, concentrations, and sample names to multiple wells with the Series feature
- Add personalized assay notes

[‡] Modes vary by instrument.

* AutoMix is covered under U.S. patent number 5,112,134. PathCheck is covered under U.S. patent numbers 5,959,738, 6,188,476, 6,320,662, 6,339,472, and 6,404,501.

[§] Available for absorbance instruments only.

Flexible and Powerful



- Apply popular linear and nonlinear regression or curve fits to recorded microplate data.
- Interpolate unknowns or data using powerful graphing options.
- Create Summary sections and enter custom formulas to summarize experimental data.
- Overlay multiple plots on a single graph, then print a report complete with raw and calculated results, graphs, and more.

FLEXIBLE CALCULATIONS AND FORMULAS

Use the extensive set of Reduction, Column, and Summary formulas, or create custom calculations to quickly analyze data.

- Perform extensive mathematical calculations within and across multiple columns of data
- Apply a single standard set across multiple plates to perform cross-plate analysis and interpolation
- Easily analyze difficult calculations like EC_{50} and Z-factors to identify potency and assay quality
- Calculate %Total Binding and %Non-specific Binding (%Control)
- Flag critical samples immediately upon reading with no manual interventions
- 4- and 5-parameter curve fit logistics and parallel line analysis (PLA)

FULL EXPERIMENTAL CONTROL

SoftMax Pro 5 Software helps organize and view experimental data in an easy, logical fashion. It offers full control to add, move, or delete sections, change font styles, and much more.

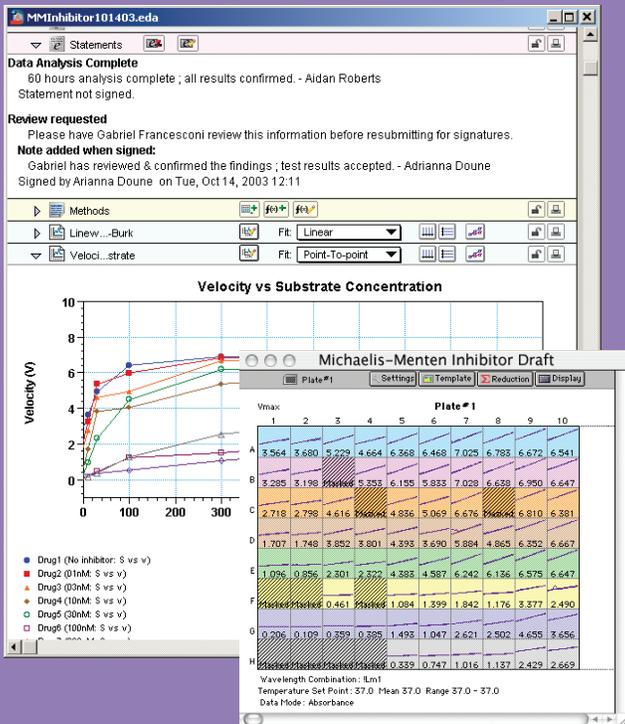
- Access curve fit parameters in formula tables
- Apply Chi-squared and F-test statistics to PLA
- Change data calculations easily using Reduction menu items
- View X/Y coordinates of any plot point
- Include multiple plates, cuvettes, graphs, analysis tables, and notes in the same data file
- Display data as OD, RFU, RLU, reduced number, grayscale, threshold, or ranged values
- Combine absorbance and fluorescence data to calculate a single result

TAILORED DATA ANALYSIS

Stop wasting time exporting data to external applications for analysis. In addition to the more than 120 predefined experiment protocols, use the straightforward yet powerful programming capabilities of SoftMax® Pro 5 software to further enhance specialized data collection and analysis results.

- Analyze data in a single plate, across multiple plates, or between different groups
- Zoom into desired wells, then adjust reduction parameters while collecting real-time data for Kinetic and Spectral Scans
- Report data using conditional formulas and apply variable weighting
- Create results summary reports

Support for FDA 21 CFR Part 11



Log On

User Name: Arianna Doune

User ID: adoune

Password: [REDACTED]

Cancel OK

Audit Trail for MMInhibitor101403.eda

Time	Date	User	User ID	Activity
8:20:34 AM	10/14/2003	User: Aidan Roberts	UserID: aroberts	User logged off
8:20:40 AM	10/14/2003	User: Guest (No Permissions)	UserID: guest	User log on attempt
8:20:52 AM	10/14/2003	User: Arianna Doune	UserID: adoune	Log on complete
8:31:28 AM	10/14/2003	User: Arianna Doune	UserID: adoune	Preferences confirmed. Instrument kind is: SpectraMax Plus 384
8:57:15 AM	10/14/2003	User: Arianna Doune	UserID: adoune	Experiment Michaelis-Menten Inhibitor: Section Statements New Statement created, Review requested Please have Gabriel Francesconi review this information before resubmitting for signatures. Not signed.
9:16:13 AM	10/14/2003	User: Arianna Doune	UserID: adoune	No user activity detected. User has been automatically logged off
12:02:39 PM	10/14/2003	User: Arianna Doune	UserID: adoune	User log on attempt
12:02:39 PM	10/14/2003	User: Arianna Doune	UserID: adoune	Log on complete
12:04:00 PM	10/14/2003	User: Arianna Doune	UserID: adoune	Document saved as M:\Customer Feedback\Customer_pda_edata files\MMInhibitor102803.eda
12:05:23 PM	10/14/2003	User: Arianna Doune	UserID: adoune	User logged off
12:05:53 PM	10/14/2003	User: Guest (No Permissions)	UserID: guest	User log on attempt
12:06:27 PM	10/14/2003	User: Guest (No Permissions)	UserID: gfrancesconi	failed password check three times. User logged off and administration file disconnected

Modify User - Craig

Permissions:

- SoftMax Pro
- SAFACS
- EBR reader settings
- EBR Note text
- Assign Plate layouts
- EBR Sample & Group Information
- Read empty plates/cuvettes
- Save data file
- ASSAY DEVELOPMENT
- ADD/DELETE GROUPS
- EBR Graphs, Summaries, and reductions
- EBR Formulas
- Lock/Unlock Sections
- Mask/UnMask wells
- Overwrite plate/cuvette data
- Save protocol files
- LAB MANAGEMENT
- Add/Modify Statements
- Change Autocore settings
- Remove Signatures
- Sign Statements

User Options:

- Force Password Change
- Active User Account
- Secondary Administrator
- Log on via Remote Command

Change the password of current user:
Change Password

Apply a predefined set of permissions:
[Dropdown]

OK Cancel

SoftMax Pro 5 GxP software contains features designed to assist a company's validation program for compliance with FDA 21 CFR Part 11 requirements including: electronic signatures, data and administrative audit trails, user management, and password aging.

POWERFUL COMPLIANCE TOOLS

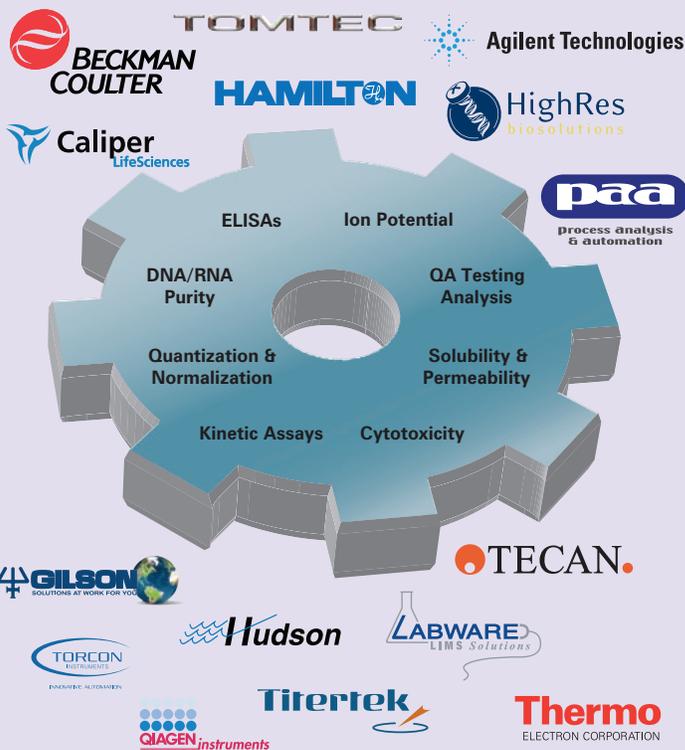
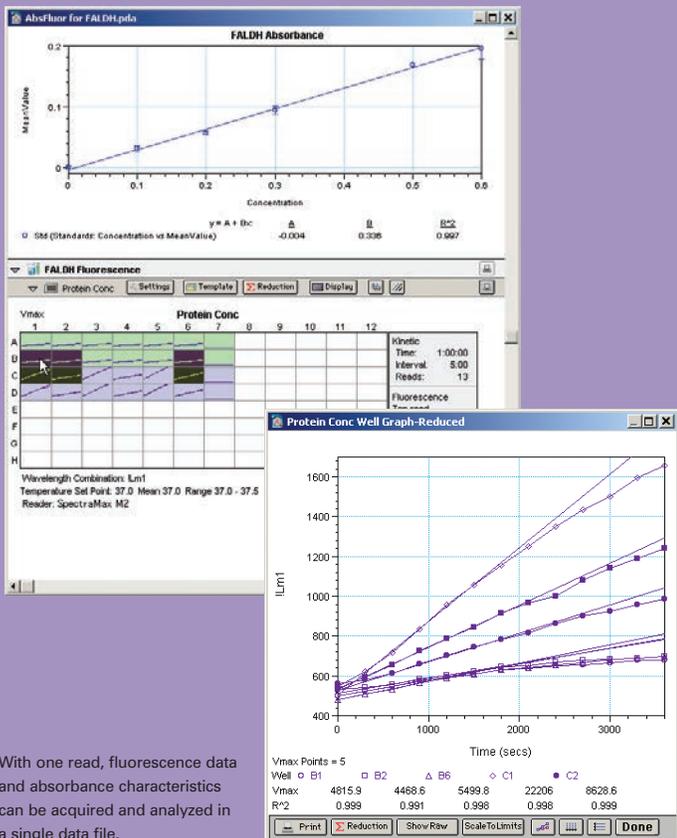
- SoftMax Pro 5 GxP Software provides FDA 21 CFR Part 11 compliance tools in conjunction with data acquisition and analysis for 23+ Molecular Devices microplate readers:
- Electronic signature support for verification, authorization, and approval
 - Powerful user management options, including multiple permission levels, templates, and more
 - Audit Trail to track and record actions
 - TCP/IP-based hosting of User Accounts files
 - AutoSave feature to save data to both local and network destinations
 - Password assignment and aging
 - Log on failure lockout and automatic idle log out
 - Centralized site-license administration

FASTER, EASIER VALIDATION

- Optional hardware and software validation tools are designed to speed and simplify the validation process by making it easier to demonstrate compliance of data collection and analysis. Available tools include:
- FDA 21 CFR Part 11 Guidance Checklist
 - Software development and quality test procedures and policies
 - Calculation, algorithm, and curve-fitting routine verification
 - SpectraTest® Validation Plates for automated SpectraMax Instrument optical verification
 - IQ/OQ/PQ validation protocols for Molecular Devices microplate readers and washers
 - In-field reader Performance Certification
 - Test Validation/Customization instructions



Widely Integrated by Molecular Devices' Automation Partners



SoftMax® Pro 5 software has been integrated by many leading robotics and LIMS partners, enabling both data analysis and instrument control in automated environments.

ROBOTICS AND LIMS INTEGRATION

- By supporting Windows-based Interprocess Messaging, SoftMax Pro 5 Software speeds and simplifies integration with leading robotics and LIMS providers and has been integrated by leading LIMS providers like LabWare and Thermo.
- Export plate data to LIMS via ASCII text or XML
 - Leverage a predefined remote command interface
 - Communicate with plate and liquid handling, plate moving, and storage robots
 - Import LIMS template information
 - Utilize abundant sample source code in Excel, MSVB, and VC++ formats

A COMPLETE PACKAGE

- For over 18 years, SoftMax Pro 5 Software has offered industry-leading support for more users than any other microplate reader software package. With a single file format that supports both Windows and Macintosh users, SoftMax Pro 5 Software is the best choice.
- Over 120 predefined assays protocols
 - Electronic and printed user documentation
 - Self-paced tutorials shorten the learning curve
 - Cross-platform file format allows easy file sharing and exchange between Windows and Macintosh users
 - Complete backwards compatibility for older SoftMax Pro Software data files

MINIMUM SYSTEM REQUIREMENTS

- Windows XP and Vista*:
- Pentium 4 1.6 GHz processor or equivalent
 - Color monitor resolution 1024 x 768
 - 512 MB of RAM
 - 90 MB of hard-disk space
 - CD-ROM drive
 - Available serial port
- Mac OS X 10.4 or later*:
- PowerPC G3 600 MHz processor
 - Color monitor resolution 800 x 600
 - 512 MB of RAM
 - 90 MB of hard-disk space
 - CD-ROM drive
 - Available 8-pin DIN serial port (ADB)
 - USB-to-Serial (ADB) adapter and software

* Windows Vista and Mac OS X 10.5 support require SoftMax Pro Software version 5.3 or greater.

SALES OFFICES

United States & Canada

Molecular Devices
Tel. +1-800-635-5577
Fax +1-408-747-3601

Brazil

Molecular Devices Brazil
Tel. +55-11-3616-6607
Fax +55-11-3871-9994

China

Molecular Devices Beijing
Tel. +86-10-6410-8669
Fax +86-10-6410-8601

Molecular Devices Shanghai

Tel. +86-21-6887-8820
Fax +86-21-6887-8890

Germany

Molecular Devices GmbH
Tel. +49-89/96-05-88-0
Fax +49-89/9-62-02-34-5

Japan

Nihon Molecular Devices Osaka
Tel. +81-6-6399-8211
Fax +81-6-6399-8212

Nihon Molecular Devices Tokyo

Tel. +81-3-5282-5261
Fax +81-3-5282-5262

South Korea

Molecular Devices Korea, LLC
Tel. +82-2-3471-9531
Fax +82-2-3471-9532

United Kingdom

Molecular Devices Ltd.
Tel. +44-118-944-8000
Fax +44-118-944-8001

www.moleculardevices.com

FOR RESEARCH ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

The trademarks used herein are the property of Molecular Devices, Inc. or their respective owners.

Specifications subject to change without notice.

©2010 Molecular Devices, Inc.
Printed in U.S.A. 6/10 #0120-1307D