



# VersaMax Microplate Reader

An affordable, tunable microplate reader for 96-well format

## BENEFITS

- Tunable for filterless flexibility
- Exceptional performance
- Affordable
- Wide application base
- Powerful data analysis
- Easy validation and testing

### Introduction

The VersaMax™ Microplate Reader from Molecular Devices bridges the gap between the affordability of filter-based readers and the flexibility of monochromator-based systems.

### Tunable, filterless flexibility

The VersaMax reader uses a grating monochromator to select the wavelength, so the exact wavelength needed can be selected for every assay. With a wavelength range from 340 to 850 nm, this is equivalent to having 510 free filters.

### Exceptional performance

Using an advanced optical and electronic design, the VersaMax reader gives the same high performance with round-bottom, flat-bottom or half-area well plates. Dual wavelength readings report the actual absorbance at each wavelength so data can be analyzed according to the user's preferences. Temperature control up to 45°C allows kinetic assays at ambient and physiological temperatures.

### Affordable

The robot-compatible VersaMax reader is cost-effective—users never have to buy filters. Furthermore, VersaMax system is reliable, which means more uptime and lower long-term cost of ownership.

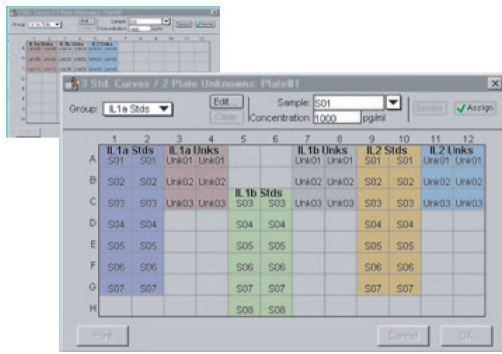
### Wide range of applications

The VersaMax reader covers a wide range of applications:

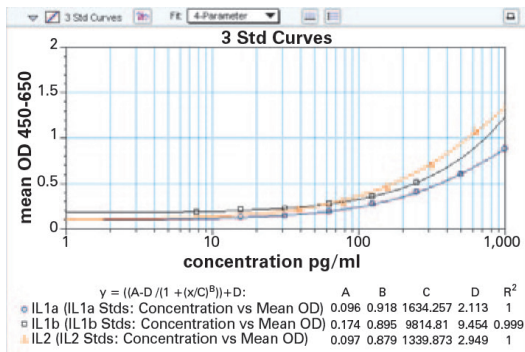
- Microbial Growth/MIC
- IC<sub>50</sub>/LD<sub>50</sub>
- Endpoint ELISAs/EIA
- Cytoproliferation/Cytotoxicity
- Colorimetric Protein
- Kinetic ELISAs/Enzyme Assays
- Bacterial Identification

### Powerful data analysis

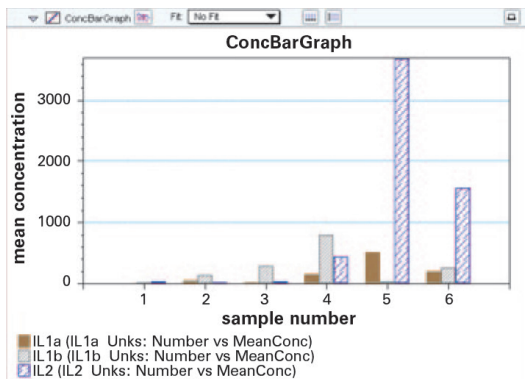
Industry leading SoftMax® Pro Data Acquisition and Analysis Software is included with VersaMax systems and provides additional flexibility for the user.



**Figure 1. Flexible template assignment.** Standards for multiple calibration curves and unknowns can be run on separate plates.



**Figure 2. Multiple calibration curves.** Multiple calibration curves can be plotted on one graph.



**Figure 3. Combined data graphing.** Results from unknowns run on different plates and different calibration curves can be plotted on one graph.

## 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.

SoftMax Pro Software can handle analysis requirements from simple endpoint assays to complex kinetic assays requiring custom calculations, meeting the needs of both basic and power users. Choose from nine different curve-fitting routines and use default data reduction, or set up custom formulas for analysis. Data can be analyzed and combined from different plates. (See Figures 1, 2 and 3.)

## Validation made easy

The SpectraTest® Validation Package tests optical performance using NIST-traceable standards. Testing can be done in the user's lab on their own schedule. Software validation, FDA 21 CFR Part 11 compliance tools and IQ/OQ/ PQ tools are also available.

### Technical specifications

#### Photometric performance

Wavelength range	340–850 nm
Wavelength selection	Monochromator, tunable in 1.0 nm increments
Wavelength bandwidth	2 nm
Wavelength accuracy	< ±2.0 nm
Photometric range	0–4.000 OD
Photometric accuracy	< ±0.006 OD ±1.0%, 0–2.0 OD
Photometric precision	< ±0.003 OD ±1.0%, 0–2.0 OD
Microplate read times	
Endpoint	12 seconds
Kinetics	9 second min. interval

#### Temperature regulation

Temperature range	Ambient +4°C to 45°C
Temperature uniformity (microplate)	±0.5°C at 37°C, well-to-well

#### General specifications

Dimensions (in.)	8.6 (L) x 22.8 (W) x 15 (H)
Dimensions (cm)	22 (L) x 58 (W) x 38 (H)
Weight	30 lbs. (13.6 kg)
Power consumption	< 250 watts
Power source	100–240 Vac, 4 A 50/60 Hz
Robot compatible	Yes

## Ordering information

Contact your Molecular Devices sales representative for configuration options.