



SpectraMax ABS and ABS Plus Microplate Readers

Small but mighty absorbance plate reader

- Affordable monochromator-based plate reader with linearity of up to 3 OD
- Works with 96- and 384-well plates as well as cuvettes (ABS Plus)
- Option to run samples as low as 2 μ L with the SpectraDrop™ Micro-Volume Microplate
- Runs various assays within the Vis/UV-Vis range such as ELISA, Bradford, BCA, microbial growth, protein quantitation and more

19,400+

Citations on Google Scholar for Molecular Devices Absorbance Microplate Readers

10,700

Citations on Google Scholar for SoftMax Pro Software

Key features

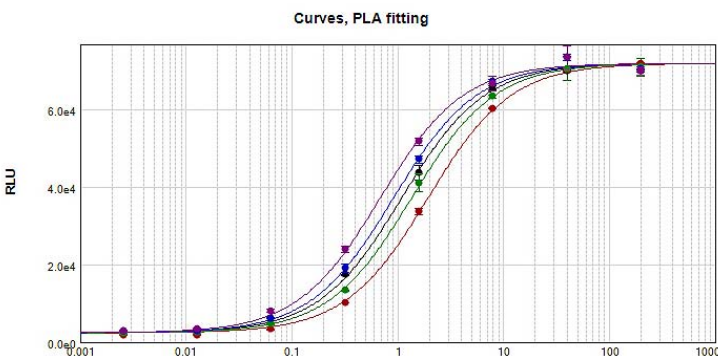
- **Convenience:**
 - Flexibility to choose your wavelength
 - Elimination of well-to-well pipetting differences with Patented PathCheck® Sensor
- **Ease-of-use:**
 - Preconfigured protocols such as ELISA and Bradford in SoftMax® Pro Software
 - Easy and fast assay setup and data analysis
- **All-around support**
 - Benefit from nearly 20,000 citations*
 - Get help with your work from our team of highly rated Ph.D. level scientists

SpectraMax ABS and ABS Plus Application Capabilities

Capabilities	SpectraMax ABS	SpectraMax ABS Plus	Wavelength	SoftMax Pro Pre-configured Protocols
ELISA	•	•	Variable	•
Bradford, BCA Protein Assay	•	•	540, 590, 595	•
NAD/NADP Conversion		•	340	•
Alkaline Phosphatase (AP) ELISA	•	•	405	•
β -galactosidase (Beta GAL)	•	•	405	•
DNA/RNA Quantification		•	260, 280	•
Horseradish Peroxidase (HRP) ELISA	•	•	492	•
MTT & XTT Cell Viability Assay	•	•	570	•
Lowry	•	•	650	•
Diphenylamine DNA (DISHE)	•	•	595	•

Data analysis made easy

- Quickly export to Microsoft Excel or utilize SoftMax Pro® Software (4 licenses included)
- 160+ preconfigured protocols (including those listed in the above table)
- Optimized analysis algorithms
- 21+ different curve fit 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