

# 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  $\mu\text{L}$  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

## Benefits

### 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:

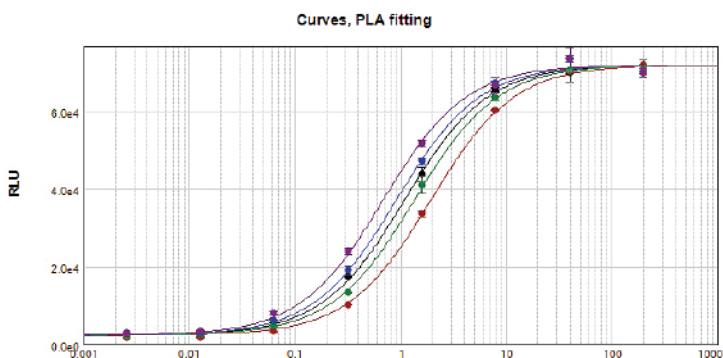
- Benefit from nearly 20,000 citations\*
- Get help with your work from our team of highly rated Ph.D. level scientists

## 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	✓
B-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](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)  
 Check our website for a current listing of worldwide distributors.

\* Source: Google Scholar search based on SpectraMax Absorbance Readers, September 17, 2018

