



# SpectraMax® QuickDrop™ Micro-Volume Spectrophotometer

Rapid, accurate DNA, RNA, and protein quantitation in a one-touch, full-spectrum micro-volume absorbance reader.

## KEY FEATURES

- Touch-screen interface with pre-configured analysis methods
- Fixed path length requires no calibration
- Micro-volume sample port has no moving parts to break or skew results regardless of viscosity
- Easily export data to a USB drive for additional analysis
- One-swipe cleaning

The SpectraMax® QuickDrop™ Micro-Volume Spectrophotometer addresses the needs of life science researchers who quantify very small amounts of DNA, RNA, oligos, and proteins. The QuickDrop Spectrophotometer allows you to set up your laboratory with a minimal investment of time, cost, and effort.

### Quantitate micro-volume samples with ease

The QuickDrop Spectrophotometer provides you with the flexibility to work with your micro-volume samples in the platform that best suits your research. The built-in sample port allows you to work with volumes as small as 0.5  $\mu$ L, preserving your precious samples, while the cuvette port lets you expand your work to include larger volume samples.

The QuickDrop Spectrophotometer is a stand-alone unit with a small footprint and does not require a direct connection to a dedicated computer. The large, high-resolution touch-screen interface offers you the ability to take advantage of pre-configured analysis methods, set up your customized experiment with ease, or export your data to a USB flash drive for additional analysis.

### Increased sensitivity without sacrificing speed

With a four second read time and no moving parts to skew the sample, or to potentially break, the QuickDrop Spectrophotometer maintains an accurate pathlength, providing you with results that are both fast and reliable, regardless of viscosity.



Quantitate micro and large volume samples in one platform using the 0.5  $\mu$ L sample port or the cuvette port.



Easy-to-use touch-screen interface allows you to easily set up your experiments or use pre-configured analysis methods in five languages.



One-swipe cleaning streamlines your workflow and allows you to quickly move from sample to sample.



Export data to your computer for additional analysis using a USB Flash Drive. No dedicated computer required.

## SpectraMax QuickDrop Micro-Volume Spectrophotometer Specifications

ITEM	DESCRIPTION	
Part Number	QUICKDROP	
Display	5.7" color, capacitive touch screen	
Light Source	Pulsed Xenon lamp	
Detector	1024 element CCD array	
	Micro-Volume Sample Port	Cuvette Port
Minimum Volume	0.5 µL	0.6 µL*
Maximum Concentration	2,500 ng/µL	10,000 ng/µL
Detection Limit (dsDNA Measurement)	1 ng/µL	0.5 ng/µL
Pathlength	0.5 mm	1 mm, 2 mm, 5 mm, and 10 mm options
Wavelength Range	190 - 1100 nm	
Wavelength Accuracy	±2 nm	
Stray Light	<0.5%T at 220 nm and 340 nm using NaNO <sub>2</sub>	
Absorbance Range	- 0.3A to 2.5A (-0.3A to 50A: 10 mm equivalent)	
Absorbance Accuracy	±0.005A or 1% of the reading, whichever is the greater at 546 nm	
Absorbance Reproducibility	±0.003A (0 to 0.5A), ±0.007A (0.5 – 1.0A)	
Output	USB port for USB Flash Drive	
Dimensions (H x W x D)	7.48 x 11.42 x 16.14 in. (190 x 290 x 410 mm)	
Weight	< 8.8 lbs. (< 4 kg)	
Display Languages	English, Simplified Chinese, French, German, Spanish	

\* varies, based on specific cuvette used



### Download Application Note

Nucleic acid quantitation and analysis using the SpectraMax QuickDrop Spectrophotometer

[moleculardevices.com/QDappnote](http://moleculardevices.com/QDappnote)



For more information on the SpectraMax QuickDrop Micro-Volume Spectrophotometer, please visit

[www.moleculardevices.com/quickdrop](http://www.moleculardevices.com/quickdrop)

### Contact Us

Phone: [+1-800-635-5577](tel:+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.