

SpectraMax Mini Multi-Mode Microplate Reader

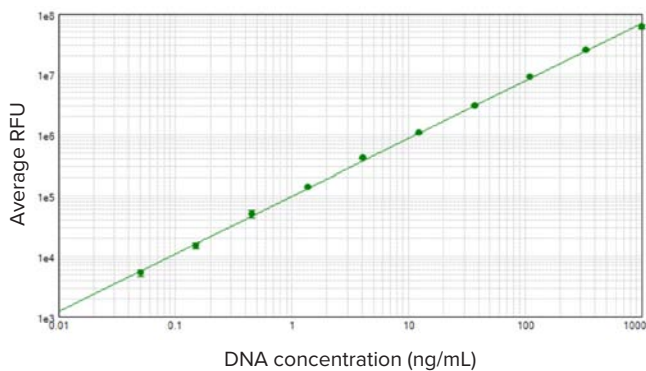
Affordable multi-mode microplate reader with best-in-class data analysis software

Perform your favorite applications including ELISA, DNA and protein quantitation, cell viability, gene expression with a user-friendly reader

With three modes of detection for full spectrum absorbance, fluorescence and luminescence, the SpectraMax® Mini reader is user upgradeable—from 2 to 3 modes, providing researchers with a versatile instrument that does not compromise on sensitivity or performance. Included SoftMax® Pro Software enables rapid data generation and analysis helping you achieve more in less time.

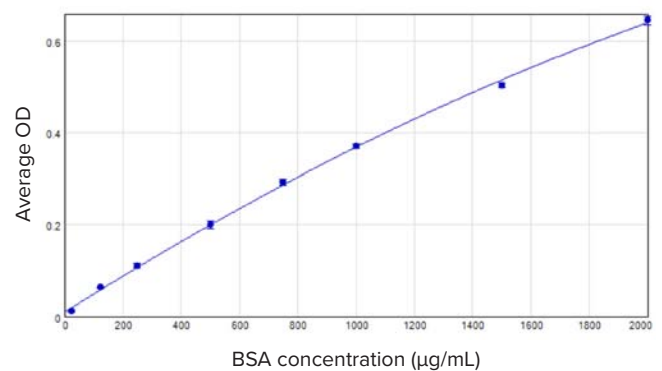
Fluorescent DNA quantitation

Determine DNA concentrations precisely and accurately over a wide dynamic range



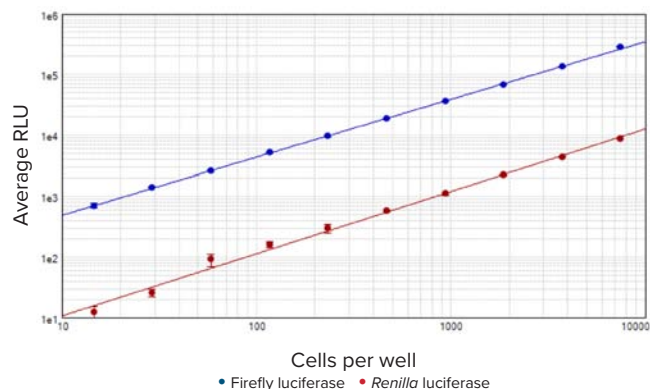
Protein quantitation

Determine protein concentrations in cell lysates and more



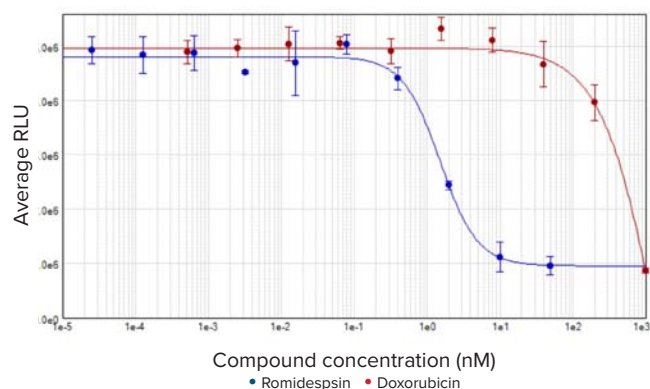
Reporter gene assays

Assess firefly and *Renilla* luciferase activity over a broad range of cell numbers and expression levels



Cell viability

Easily measure cell viability with a luminescence readout



General specifications

Dimensions (in.)	15.75 (L) x 12.60 (W) x 13.78 (H)
Dimensions (cm)	40 (L) x 32 (W) x 35 (H)
Weight	≤ 40 lbs (18 kg)
Power requirement	External 24V DC power supply compatible with 100–240V AC @50–60 Hz 160W maximum consumption
Robotic compatible	Yes

General performance

Plate formats	6, 12, 24, 48, 96, 384-well plates ANSI/SLAS
Light source	Two high-energy xenon flash lamps
Reading capabilities	Microplates, cuvettes (via adapter)
Detectors	Ultra-cooled Photomultiplier Tube and Silicon photodiode
Shaking	Linear, orbital, and double orbital
Temperature control	4°C above ambient to 45°C
Temperature uniformity	±1°C at 37°C
Temperature accuracy	±2°C at 37°C
Spectral scanning	Abs
Endpoint reading	Abs, FI, Lum
Kinetic reading	Abs, FI, Lum
Well scanning	Over 20 by 20

Absorbance photometric performance

Wavelength range	200–1000 nm
Wavelength selection	1.0 nm increments
Wavelength bandwidth	≤ 5 nm full width half maximum
Wavelength accuracy	± 2 nm across wavelength range
Wavelength precision (repeatability)	± 0.2 nm
Photometric range	0–4.0 OD
Photometric resolution	0.001 OD
Photometric accuracy, linearity	< ±0.006 OD ±1.0%, 0–3 OD
Photometric precision (repeatability)	< ±0.003 OD ±1.0%, 0–3 OD
Stray light	≤ 0.05% at 230 nm

Fluorescence intensity performance

Wavelength range	250–850 nm
Dynamic range	> 6 decades
Sensitivity (fluorescein)	1 pM

Luminescence performance

Wavelength range	300–650 nm
Dynamic range	> 7 decades
Sensitivity (ATP)	2 pM
Cross-talk	<0.1% (white 96-well half-area microplate)

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
 United Kingdom +44.118.944.8000
 Europe* 00800.665.32860
 China +86.4008203586
 Taiwan/Hong Kong +886.2.2656.7585
 Japan +81.3.6362.9109
 South Korea +82.2.3471.9531
 India +91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom