



Absorbance

SpectraMax Detection Cartridge

Absorbance SpectraMax® Detection Cartridges for the SpectraMax® Paradigm® System combine scanning and a broad spectrum wavelength range necessary to address a variety of nucleic acids, proteins, ELISAs and immunoassays found in the laboratory.

Absorbance SpectraMax Detection Cartridge

Description

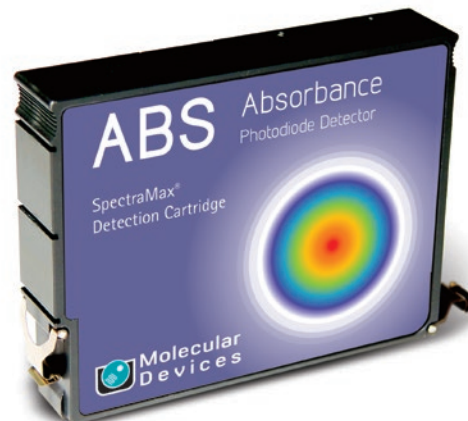
Compatible system	SpectraMax Paradigm Reader
Part Number	0200-7000
Mounting	Top Read (Photodiode Detector) and Bottom Read (Monochromatic Illumination)
Cartridge Size	Top - 1 Position Bottom - 2 Position
Light Source	Flash Lamp

Technical specifications

Wavelength Range	230-1000 nm
Wavelength Selection	Monochromator (1.0 nm increments)
Wavelength Bandwidth	4.0 nm
Wavelength Accuracy	± 1 nm
Wavelength Repeatability	± 0.5 nm
Photometric Range	0 - 3.5 OD
Photometric Resolution	0.0001 OD
Photometric Accuracy	$< \pm (2\% + 0.010 \text{ OD})$, 0-2 OD @ 405 nm
Photometric Precision	$< \pm (1\% + 0.005 \text{ OD})$, 0-2 OD @ 405 nm

Key applications

- Nucleic Acid Quantitation
- Protein Quantitation
- ELISA
- Immunoassay
- Proliferation/Viability
- Low Volume Applications



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	Japan (Osaka)	+81-6-7174-8831
Brazil	+55-11-3616-6607	Japan (Tokyo)	+81-3-6362-5260
China (Beijing)	+86-10-6410-8669	South Korea	+82-2-3471-9531
China (Shanghai)	+86-21-3372-1088	United Kingdom	+44-118-944-8000
Germany	00800-665-32860		

