

SpectraMax Paradigm

Absorbance Detection Cartridge



The **SpectraMax® Paradigm® Absorbance Detection Cartridge** combines scanning and a broad spectrum wavelength range necessary to address a variety of nucleic acids, proteins, ELISAs and immunoassays found in the laboratory.

SpectraMax Paradigm Multi-Mode Detection Platform

Absorbance Detection Cartridge

DESCRIPTION

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

TYPICAL PERFORMANCE

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

TYPICAL APPLICATIONS

- Nucleic Acid Quantitation
- Protein Quantitation
- ELISA
- Immunoassay
- Proliferation/Viability



Contact Us

Phone: +1-800-635-5577
Web: www.moleculardevices.com
Email: info@moldev.com
Check our web site for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1-800-635-5577
Brazil +55-11-3616-6607
China (Beijing) +86-10-6410-8669
China (Shanghai) +86-21-6887-8820
Germany +49-89/96-05-88-0

Japan (Osaka) +81-6-6399-8211
Japan (Tokyo) +81-3-5282-5261
South Korea +82-2-3471-9531
United Kingdom +44-118-944-8000