



xMark™ Microplate Absorbance Spectrophotometer

Extending the Limits of Microplate Reading With Remarkable Flexibility

With its unique imaging capabilities, monochromator design, and spectral scanning feature, the xMark microplate absorbance spectrophotometer finds the best wavelength for any photometric application, and supports assay development in a wide variety of formats.



| Feature | Benefit |
|--|--|
| Unique monochromator design with wide spectral range of 200–1,000 nm and 1 nm step size | <ul style="list-style-type: none"> Detection, analysis, and quantitation of substances with absorbance spectra in the UV range (DNA) to the infrared range (water) |
| Capability to scan standard 6- to 1,536-well microplates and custom plates | <ul style="list-style-type: none"> High throughput with sample volumes below 5 µl No restrictions on sample type, volume, or throughput Wide range of applications, from determining HLA antigens serologically to testing wine quality |
| Unique well and plate imaging | <ul style="list-style-type: none"> Intensity maps for analysis of absorbance variations in a single well or entire plate Enhanced detection sensitivity |
| Built-in incubator* with continuous temperature display using Microplate Manager® software | <ul style="list-style-type: none"> Thermal regulation of assay reactions Improved sensitivity and activity of assays Incubation of plates prior to detection |
| Built-in shaker* with variable mixing speed and linear or orbital mixing mode | <ul style="list-style-type: none"> Accurate absorbance readout for entire well volume Minimum variability of assays for different sample types and liquid- or solid-state reactions Assay preparation and repeated mixing |
| Single and dual wavelength detection | <ul style="list-style-type: none"> Detection of reaction products with high specificity while eliminating background and noise |
| Motorized tray for plate loading and retrieval | <ul style="list-style-type: none"> Minimized environmental exposure |
| Fast plate reading — 9 sec at single wavelength detection for a 96-well plate | <ul style="list-style-type: none"> Precision in kinetic assays and rapidly developing colorimetric reactions |
| Microplate Manager software controlled (PC or Mac) via USB2 port | <ul style="list-style-type: none"> Protocol customization and data reduction options Comprehensive data analysis |
| Reports with Checkmark™ reader performance validation kit | <ul style="list-style-type: none"> Accommodation of various validation protocols, including installation qualification and operational qualification (IQ/OQ) |

* Incubator and shaker can be used independently from other instrument functions.



Specifications

| Optical | |
|-------------------------|---|
| Wavelength | |
| Range | 200–1,000 nm |
| Accuracy, minimum | ±2.0 nm |
| Repeatability | ±2.0 nm |
| Photometric methods | Single and dual wavelength |
| Monochromator step size | 1 nm |
| Bandwidth | 5 nm |
| Imaging resolution | 0.5, 0.75, 1.0 mm |
| Performance | |
| Indication range | 0–4.0 OD |
| Resolution | 0.001 OD |
| Linearity | ≤2.0%, 0–3.0 OD at 405 nm |
| Accuracy | ±1.0% or 0.015, 0–3.0 OD at 405 nm |
| Reproducibility | 96-well plate; ≤1.0% or 0.005, 0–3.0 OD at 405 nm |
| Stability and drift | ≤0.010 at OD = 1 at 490 nm |
| Read time | 8 sec/96-well/single wavelength; 15 sec/384-well/single wavelength; 35 sec/1,536-well/single wavelength |
| Physical | |
| Compatible plates | 6- to 1,536-well and custom microplates; maximum plate height: 21 mm |
| Incubator set point | Temperature range: 25–45°C; accuracy: ±0.5°C; uniformity: 0.5°C |
| Well-to-well uniformity | ≤2.0% or 0.005, 0 to 3.0 OD at 405 nm |
| Plate shaking | Linear or orbital plate-shaking capability |
| Computer interface | USB2 |
| Dimensions (W x D x H) | 54.4 x 44.0 x 21.9 cm (21.4 x 17.3 x 8.6") |
| Weight | 27.5 kg (61 lb) |

Ordering Information

| Catalog # | Description |
|-----------|--|
| 168-1150 | xMark Microplate Absorbance Spectrophotometer , PC or Mac, includes incubator, Microplate Manager software, USB2 and power cables |

Mac is a trademark of Apple Inc.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400 **Canada** 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000