



iMark™ Microplate Absorbance Reader

On the Mark for Cost-Effective and Reliable Absorbance Detection

The iMark microplate absorbance reader is an economical, high-performance solution for a wide range of photometric applications. The iMark reader represents the next generation of Bio-Rad's popular Model 550 and Model 680 series microplate readers. For over two decades, Bio-Rad absorbance microplate readers have hit the highest marks for reliability and efficiency in thousands of laboratories around the world.



Feature	Benefit
Onboard software, featuring an intuitive user interface, preset or customized end-point and kinetic analysis protocols, powerful curve-fitting routines, and statistical analysis	<ul style="list-style-type: none"> Automatic control of detection, data formats, and analyses Data storage and reports for most complex photometric applications Easy configuration, storage, and application of customized protocols
Test protocols and reporting for transmissible spongiform encephalopathies (TSEs), such as BSE (bovine)	<ul style="list-style-type: none"> Minimum intervention during testing of food-related samples
Integrated thermal printer	<ul style="list-style-type: none"> Instant documentation of real-time data
Automatic calibration before each reading	<ul style="list-style-type: none"> Accurate and reliable measurements
Built-in, variable-speed plate-shaking capability	<ul style="list-style-type: none"> Accurate absorbance readout for entire well volume Minimum variability of assays for different sample and reaction types
Ability to read flat-, U-, or V-bottom microplates or 8- or 12-well strip plates	<ul style="list-style-type: none"> Optimum assay and reaction conditions and sample recovery
Single and dual wavelength detection	<ul style="list-style-type: none"> Detection of reaction products with high specificity, while eliminating background and noise
Motorized door for plate loading and retrieval	<ul style="list-style-type: none"> Minimum environmental exposure
Multilingual interface, LCD display, and reports (English, Japanese, Chinese, and Russian)	<ul style="list-style-type: none"> Accommodation of localized regulatory requirements and communication standards
Fast plate reading — 6 sec at single wavelength detection	<ul style="list-style-type: none"> Rapid detection for precision in kinetic assays and rapidly developing colorimetric reactions
Fast or step modes	<ul style="list-style-type: none"> Fast reading or high-sensitivity detection
Self-diagnostic capabilities to detect lamp burnout at startup	<ul style="list-style-type: none"> Proactive and prompt maintenance
USB2 port for optional external computer control (with Microplate Manager® software, PC or Mac)	<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)

BIO-RAD



iMark Microplate Absorbance Reader

Specifications

Wavelength range	400–750 nm
Photometric methods	Single or dual wavelength
Photometric range	0.0–3.5 OD
Bandwidth	10 nm
Linearity	≤1.0% from 0.0–2.0 OD; ≤2.0% from 0.0–3.0 OD
Accuracy	±1.0% or 0.010 from 0.000–3.000 OD at 490 nm
Reproducibility	1.0% or 0.005 OD from 0.0–2.0 OD; 1.5% from 2.0–3.0 OD
Resolution	0.001 OD
Light source	Tungsten halogen lamp (20 W), 3,000 hr average lifetime
Photodetectors	Silicon photodiodes: 8 measurement, 1 reference
Filters	8-position filter disk
Read time	Fast mode: 6 sec at single wavelength, 10 sec at dual wavelength Step mode: 15 sec at single wavelength, 25 sec at dual wavelength
Plate shaking	3 speeds: low, mid, high; duration: 0–999 sec
Plate types	96-well microplate; maximum plate height: 16 mm
Warm-up time	3 min
Interchannel variation	≤1.0% or 0.005 from 0 to 3.0 OD
Stability and drift (at 490 nm)	≤0.010 OD at OD = 1 at 490 nm single wavelength
Data output	Onboard graphical thermal printer and USB2 interface with PC or Mac data stations
Data storage	Calendar/clock function; 64 assay protocols
Multilanguage support	4 language ROM capacity; LCD indication supported; printout report supported
Dimensions (W x D x H)	34.6 x 37.7 x 16.4 cm (13.6 x 14.8 x 6.5")
Weight	5.5 kg (12 lb)

Ordering Information

Catalog #	Description
168-1130	iMark Microplate Absorbance Reader , 100/240 V, includes 4 filters (415, 450, 490, 655 nm), plate shaker, onboard software and thermal printer, one roll printer paper, USB2 and power cables, instructions
168-1135	iMark Microplate Absorbance Reader With Microplate Manager Software

Mac is a trademark of Apple Inc.


BIO-RAD

**Bio-Rad
Laboratories, Inc.**



Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670
The Netherlands 0318 540666 **New Zealand** 64 9 415 2280 **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 3170 **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 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000