



# PR 4100 Absorbance Microplate Reader

## The Next Generation Reader

The PR 4100 is the next generation 8-channel absorbance microplate reader by Bio-Rad. Along with its compact and innovative design, it provides accurate, reproducible and fast measurements. Controlled by a Notebook PC with Windows 7-based software, the system is validated with the complete menu of Bio-Rad EIA tests.

### Improved Reliability

- Latest LED Technology for long life and maintenance free operation
- 8-channel absorbance microplate reader for fast measurements
- Comprehensive QC function with multicheck plate

### Innovative Design

- Compact footprint
- Standalone reader operated through a Notebook PC
- Windows® 7-based software

### Comprehensive Data Analysis Software

- Validated protocols according to Bio-Rad quality assurance process
- Power Data Analysis with full qualitative and quantitative EIA analysis
- Full traceability

### New LIS Connectivity

- Transfer to Laboratory Information Service for all Bio-Rad tests
- ASTM standard protocol

**BIO-RAD**

## PR 4100 Absorbance Microplate Reader

### Specifications

Reading head . . . . . 8 measurements, 1 reference channel  
 Light source . . . . . LED, Self-calibrating  
 Measurement range . . . . . 0-4.0 OD  
 Wavelength range . . . . . 400-750 nm  
 Wavelength selection . . . . . Interference filters (wheel of 8 filters)  
 . . . . . Supplied: 405, 450, 492, 550 and 620 nm  
 . . . . . Bandwidth: 10 nm at 50% transmission  
 Measurement time:  
 Single wavelength . . . . . < 15 sec  
 Dual wavelength . . . . . < 20 sec  
 Accuracy (450/492 nm):  
 0.000 – 2.000 OD . . . . . Better than ± (1.0% + 0.010 OD)  
 2.000 – 3.000 OD . . . . . Better than ± (1.5% + 0.010 OD)  
 Precision (450/492 nm):  
 0.000 – 2.000 OD . . . . . Better than ± 1.0%  
 2.000 – 3.000 OD . . . . . Better than ± 1.5%  
 Linearity (450/492nm):  
 0.000 – 2.000 OD . . . . . Better than +/- 1.0 %  
 2.000 – 3.000 OD . . . . . Better than +/- 1.5 %  
 Resolution . . . . . 0.0001 OD  
 Shaking . . . . . Linear shaking in 4 different modes

### Computer & Software

Operating system . . . . . Windows® 7  
 Software . . . . . Magellan Tracker

### Physical Data

Size in Inches (H x W x D) . . . . . 5.3 x 13.7 x 7.4  
 Weight . . . . . 5.7 lbs

### Electrical

Main power input . . . . . Basic instrument with AC adapter  
 External power supply . . . . . 100-240V AC: 50/60 Hz (Auto sensing), max 1.2 A  
 Power consumption:  
 Standby mode . . . . . Approx 12 VA  
 Operational mode . . . . . Max 30 VA

### Software Language

English

### Optional

Wide range of filters  
 Calibration test plate with software (Multicheck plate)

## PR 4100 is Part of a Complete System for ELISA Tests



PR 4100 Absorbance  
Microplate Reader



PW41 Microplate Washer



IPS Microplate Incubator

## Ordering Information

Catalog No.	Description	Quantity
45591587	PR 4100 Absorbance Microplate Reader (includes AC Adaptor)	1 unit



Bio-Rad  
Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics)

Clinical  
Diagnostics Group

Website [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) U.S. 1-800-224-6723 Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3237-9400  
 Canada 1-514-334-4372 China 86-21-61698500 Czech Republic 420-241-430-532 Denmark +45-4452-1000 Finland 358-9-804-22-00 France 33-1-47-95-60-00  
 Germany +49-(0)89-318-840 Greece 30-210-7774396 Hong Kong 852-2789-3300 Hungary +36-1-459-6100 India 1800-180-1224 Israel 972-3-9636050 Italy +39-02-216091  
 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico +52 (55) 5488-7670 The Netherlands +31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30  
 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia 7-495-721-14-04 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200  
 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44-(0)20-8328-2000