



# PR 3100 TSC Microplate Reader

## Consolidated testing and reporting that meets your needs

The PR 3100 TSC Microplate Reader is a stand-alone, computer-driven reader which measures the absorbance of specimens in 96-well plates. On-board data reduction generates accurate cutoff calculations after analyzing data from endpoint measurements. The wavelength range of 402 to 620 nm ensures compatibility for the broad, validated, testing menu on the reader. The ability to connect a barcode reader, printer and USB storage device for data export increases the functionality of this microplate reader.

### Flexibility

- Data export to Excel or LIS
- Single and dual wavelength reading
- Automated sample ID

### Consolidation

- Broad testing capacity
- Comprehensive data reduction

### Compact

- On-board computer
- ~ 4000 plate result storage



**PR 3100 TSC Microplate Reader**

**Instrument Specifications**

**Electrical**

Line Voltage .....	100-120 and 220-240 V ± 10%
Frequency .....	50/60 Hz
Power Consumption .....	110 VA

**Physical**

Dimensions .....	13.2" wide x 12.6" diameter x 6.7" high (33.5 cm x 32 cm x 17 cm)
Weight .....	13.2 lbs (6 kg)

**Environmental**

Temperature .....	59-95°F (15-35°C)
Relative Humidity .....	20-90%

**Data Management**

Operating System .....	Windows CE
Software .....	Magellan CE v5.2
Storage 1 .....	256 MB (APF USB memory stick)
Storage 2 .....	256 MB (Export USB memory stick) – ASTM & ASCII
Display .....	Colored Touchscreen – 640 x 480 pixels

**Optical**

Light Source .....	Halogen Lamp 20W
Measurement Range .....	0-4.0 Abs
Wavelength Range .....	340-750 nm
Wavelength Selection .....	Interference filters (slide up to 6 filters)
Filters supplied .....	402, 450, 492, 550, 620 nm
Reading Speed .....	6 sec (single), 8 sec (dual)
Resolution .....	0.001 Abs

<b>Accuracy</b>	492 nm	0.000-2.000 Abs: Better than ± 1.0% ± 0.010 Abs
		2.000-3.000 Abs: Better than ± 1.5% ± 0.010 Abs
<b>Precision</b>	492 nm	0.000-2.000 Abs: Better than ± 0.5% ± 0.005 Abs
		2.000-3.000 Abs: Better than ± 1.0% ± 0.005 Abs
<b>Linearity</b>	340-399 nm	0.000-2.000 Abs: Better than ±2%
	400-750 nm	0.000-2.000 Abs: Better than ± 1%

**Ordering Information**

Catalog No.	Description	
92000B	PR 3100 TSC Absorbance Microplate Reader (includes Magellan v5.2 software and User Manual) .....	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. 1-800-2BIO-RAD 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 86-21-64260808 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 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216091  
Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico 52(55)5200-0520 The Netherlands +31-318-540686 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 65-6415-3188 South Africa 27-11-442-85-08 Spain 34-91-590-5200  
Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom +44-(0)20-8328-2000