



SmartSpec™ Plus spectrophotometer



Simply Brilliant.

Performance, stability, and functionality at an affordable price.

BIO-RAD

The SmartSpec Plus

spectrophotometer



Features

- Compact, space-saving design
- Built-in printer records results immediately
- Easy-to-use, menu-driven operation
- Xenon flash lamp extends lamp life and reduces maintenance costs
- Capable of spectral scans from 200 to 800 nm
- Multilanguage interface — English, French, German, Italian, Spanish, and Japanese display

Accessories

Quartz Cuvettes

Bio-Rad provides a selection of quartz cuvettes for your nucleic acid assay needs. All quartz cuvettes are heat fused and annealed to relieve strain. Bio-Rad quartz cuvettes will resist most solvents and mild acidic and basic solutions. Choose from among four cuvette sizes.

Protein Assay Cuvettes

Bio-Rad maintains an inexpensive line of disposable cuvettes for protein assays. Both the 3.5 ml and 1.5 ml cuvettes are supplied in packages of 100 and are compatible with the SmartSpec Plus.

Colorimetric Protein Assays

Bio-Rad offers three colorimetric assays for protein quantitation: the Bio-Rad protein assay for general use, the *DC*[™] protein assay for samples solubilized in detergent, and the *RC DC*[™] protein assay for samples in the presence of both reducing agents and detergents.

Performance Validation Package

The SmartSpec Plus may be purchased with an optional performance validation package that allows routine checks of critical performance specifications. This package includes a NIST-traceable reference cuvette for checking wavelength and photometric accuracy, and a certificate of analysis.

Coming Soon!

trUView[™] Cuvettes

Bio-Rad will be offering low-volume (50 µl), disposable plastic cuvettes suitable for assays measured between 220 and 1,100 nm. A single-use alternative to quartz cuvettes, trUView cuvettes are designed for the quantitation of small-volume nucleic acid samples. trUView cuvettes are provided in packs of 50 or 100 individually wrapped cuvettes.

Specifications

Light source	Xenon lamp
Detector	Photodiode array
Display	2 x 24 alphanumeric backlit LCD
Printer	Built-in alphanumeric/graphic printer
Interface	RS-232 serial port
Wavelength range	200–800 nm
Wavelength accuracy (at 18–27°C)	<±1.0 nm from 200 to 250, 300 to 800 nm <±0.5 nm from 250 to 300 nm
Spectral bandwidth	<5 nm
Absorption accuracy	±0.01 at 0.5 AU; ±0.02 at 1.0 AU
Absorption repeatability	±0.005 at 0.5 AU
Stray light	<0.2% at 230 nm using NaI standard
Operating temperature	15–35°C
Dimensions	33 x 26 x 14 cm (W x D x H)
Weight	~3.8 kg
Power requirements	90–260 VAC, 47–63 Hz
Regulatory	CE, TÜV

Ordering Information

Catalog #	Description
-----------	-------------

Spectrophotometer Configurations

170-2525	SmartSpec Plus Spectrophotometer
170-2520	SmartSpec Plus Spectrophotometer, includes validation package

Accessories

170-2509	Performance Validation Package, includes NIST-traceable cuvette, certificate of analysis
170-2502	Standard Cuvette, 3.5 ml, quartz
170-2503	Semimicrovolume Cuvette, 1.4 ml, quartz
170-2504	Microvolume 700 µl Cuvette, quartz
170-2505	Submicrovolume 100 µl Cuvette, quartz
223-9950	Standard Disposable Polystyrene Cuvettes, 3.5 ml, 100
223-9955	Semimicro Disposable Polystyrene Cuvettes, 1.5 ml, 100
170-2506	Printer Paper, 5-pack
170-2507	Spring, cuvette holder

Protein Assay Kits

500-0001	Bio-Rad Protein Assay Kit I, bovine γ-globulin standard, based on the method of Bradford
500-0002	Bio-Rad Protein Assay Kit II, bovine serum albumin standard, based on the method of Bradford
500-0111	<i>DC</i> Protein Assay Kit I, bovine γ-globulin standard
500-0112	<i>DC</i> Protein Assay Kit II, bovine serum albumin standard
500-0121	<i>RC DC</i> Protein Assay Kit I, includes <i>RC</i> reagents package, <i>DC</i> protein assay reagents package, bovine γ-globulin standard
500-0122	<i>RC DC</i> Protein Assay Kit II, includes <i>RC</i> reagents package, <i>DC</i> protein assay reagents package, bovine serum albumin standard

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191
Canada (905) 712-2771 Czech Republic + 420 2 41 43 05 32 China (86-21) 63052255 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127
Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54
The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 8126 672 Portugal 351-21-472-7700
Russia 7 095 721 1404 Singapore 65-62729877 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700
Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000