



Elite™ Microplate System

Automated Microplate Workstation

A solution for the high volume ELISA microplate laboratory



Elite™ Microplate System

Sample Processing

Deck Processing 54 tracks
 Allows Combinations of Max. 30 tube carriers and 1 carrier of 36 plates
 Tube Carrier Capacity 24 or 32 tubes
 Tube Capacity Up to 960/run
 Useable Diameter Up to 18 mm
 Useable Height Up to 120 mm
 Container Mix Yes
 Minimum Sample Dilution (Direct) 1 to 101
 Carry-Over None

Robotic Positioning 0.1 mm position X-Y-Z resolution

Disposable Tips

On-Board Capacity 750 tips, 300 µL and 190 tips, 1000 µL
 Unload Option Yes
 Reload Option Yes
 Waste Capacity > 1000 tips
 Tip Detection Yes (mechanical)

Pipetting Specification

Standard 300 µL
 High 1000 µL

Pipet Grade/Size	Trueness (%)	Precision (%)
Standard 10 µL	5.0	2.25
Standard 50 µL	2.0	0.75
Standard 200 µL	1.0	0.75
High 10 µL	7.5	3.5
High 50 µL	2.0	0.75
High 200 µL	1.0	0.75

Patient Samples (100 µL/well) < 13 minutes/plate
 Reagent (100 µL/well) < 4 minutes/plate

Liquid Detection Capacitive/Pressure LLD Monitoring

Autoload Bar Code Scanner

During Carrier Loading the Following
 Bar codes are Read Plates, carrier type, sample tubes,
 control tubes, reagent container
 Positive Samples/Reagent ID Yes
 Reads ISBT Standard, Code 128, Code 39, Codabar

Reader

Capacity 1 plate
 Reading Head 8 channels
 Number of Ref. Channels 1
 Filters Supplied as Standard w/ Instrument 5 (340, 405, 450, 492 & 620 nm)
 Linear Measurement Range, 400-750 nm 0.00-2.5 OD
 Indication Range 0.0-2.999 OD

Data Processing

User Interface Software under Windows® XP
 Computer Supplied
 Communication On-line (uni-directional) ASTM or ASCII format

iSWAP

Plate Format Plate height < 43 mm
 Transport Area Whole instrument work area

Plate Processing

16/20 Configuration 1 washer/dispenser,
 16 reagent positions, 20 incubator slots
 24/20 Configuration 2 washer/dispensers,
 24 reagent positions, 20 incubator slots

Incubation

Slots Individually Monitored 20
 Temperature 5 ambient, 15 ambient to 45°C

Reagent Processing

Volume of Syringe 20 mL
 Dispense Volume Per Well 20-200 µL
 In Steps of 5 µL
 Container Capacity 100 mL (approx.)
 Reagent Identification By bar code

Washer

Wash Volume 50-400 µL (Step 10 µL)
 Residual Volume After Washing Mean < 10 µL/well, typically 5 µL/well
 Number of Wash Cycles 1-9 cycles
 Wash Modes Strip wash, Plate wash
 Soak Time 20-300 seconds
 Plate Wash 0-60 seconds
 Strip Wash Continuous wash
 Wash Options Bottom wash or sweep
 Number of Wash Containers 3 per washer
 Number of Rinse Containers 1 per washer
 Wash/Rinse Container Capacity 3 liters
 Waste Container Capacity 10 liters (exchangeable during operation)
 External Water Separator Capacity 4 liters

Electrical

Line Voltage 100-260V
 Frequency 47-63 Hz
 Input Power Max. 0.9 kVA

Physical Dimensions

Height: 1670 mm (66 in.)
 Length: With 16/20 Approx. 2930 mm (115 in.)
 With 24/20 Approx. 3140 mm (124 in.)
 Depth: Without Loading Trays 806 mm (32 in.)
 With Loading Trays and Carriers 1210 mm (48 in.)
 Weight: With 16/20 Approx. 700 kg (1543 lb.)
 With 24/20 Approx. 750 kg (1653 lb.)

NOTE: Specifications are subject to change without prior notice. Elite is a trademark of Bio-Rad Laboratories. Windows is a registered trademark of Microsoft Corporation

Ordering Information

Catalog No.	Description
26173	Elite™ Microplate System 1 system



**Bio-Rad
 Laboratories**

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

*Clinical
 Diagnostics Group*

Website 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** 36-1-455-8800 **India** 91-124-4029300 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-6902 **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** 65-6415-3188 **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