



EVOLIS™ System

One system, many opportunities





One System...

Today, large laboratories processing multiple tests need to do more with less in order to stay competitive. Labor represents a large cost, so being able to organize staff efficiently is an advantage. The EVOLIS™ System provides that advantage: the operator loads samples and consumables onto each EVOLIS™ System, then walks away to work on other tasks while specimens are processed.



Trusted

- Proven solution
- Uncompromised safety



Flexible

- Adaptable to individual laboratory requirements
- Variable configuration with multiple EVOLIS™ Systems



Consolidation

- Single platform for all testing
- Comprehensive range of assays



Efficient

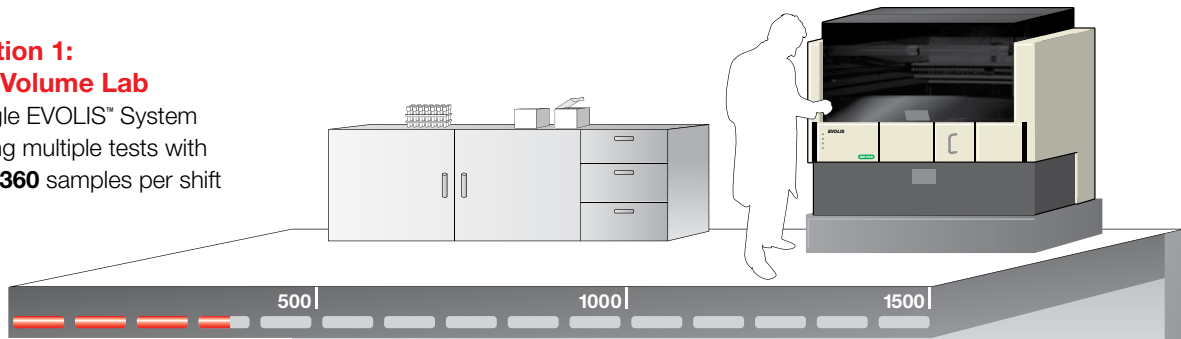
- Intuitive processes
- Full automation

Many Opportunities

With the EVOLIS™ System, you can always find the right solution

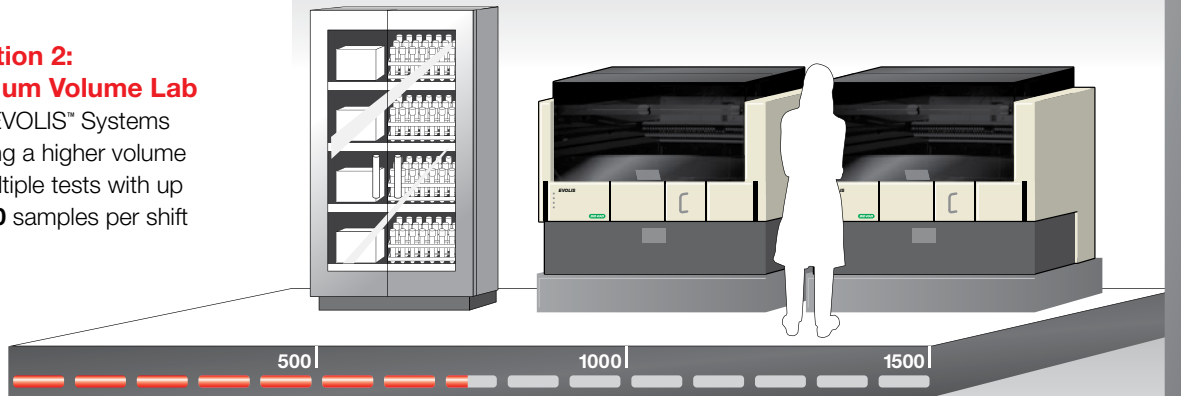
Solution 1: Low Volume Lab

A single EVOLIS™ System running multiple tests with up to **360** samples per shift



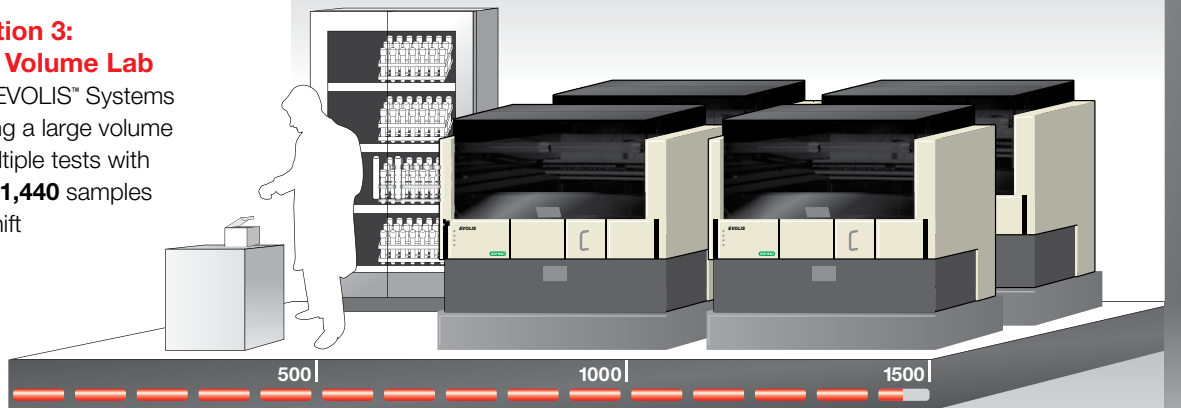
Solution 2: Medium Volume Lab

Two EVOLIS™ Systems running a higher volume of multiple tests with up to **720** samples per shift



Solution 3: High Volume Lab

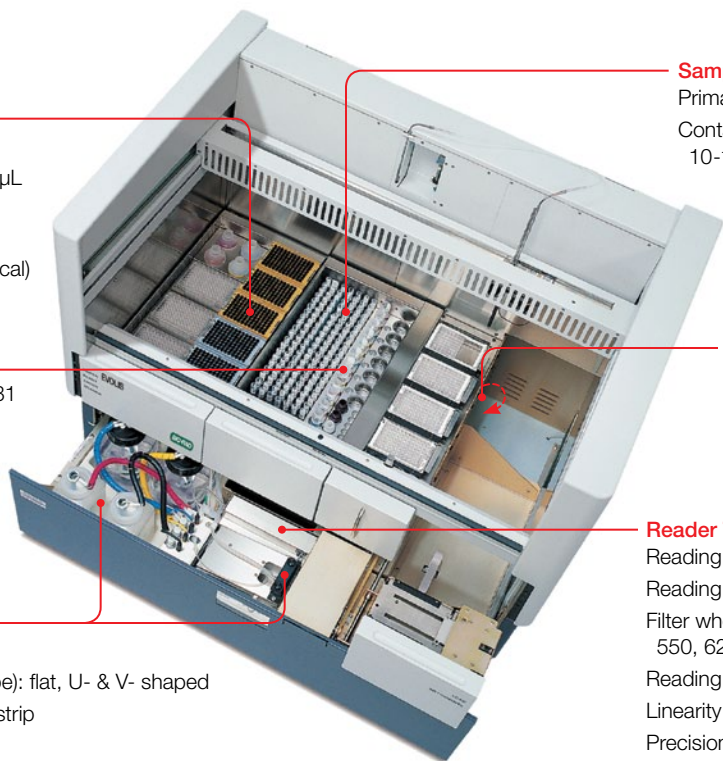
Four EVOLIS™ Systems running a large volume of multiple tests with up to **1,440** samples per shift



Expandable up to 6 EVOLIS™ Systems

EVOLIS™ System

Specifications



Disposable Tips

Carbon tips: 300-1000 µL
 Volume (useable): 10-1000 µL
 Liquid level & clot detection: capacitive
 Tip detection: yes (mechanical)

Reagent Processing

Calibrator & Control capacity: 31
 Reagent container: Bio-Rad reagent bottles
 Reagent bottle volume (mL): 3, 8, 15, 30, 60, 125

Washer Tool

Manifold: 8 channels
 Plate type (bottom shape): flat, U- & V- shaped
 Wash mode: plate and strip

Sample Processing

Primary sample capacity: 180
 Container (diameter & height): 10-16 mm/50-100 mm

Incubation Tower

Loading slots: 4 x ambient T°
 Chambers (independent): 4 x RT + 5-50°C

Reader Tool

Reading head: 8 channels
 Reading range: 0-3.5 O.D.
 Filter wheel: 8 (equipped 405, 450, 492, 550, 620, 690 nm)
 Reading method: (beam) single/double
 Linearity (0-2.0 O.D.): 1%
 Precision (0-2.0 O.D.): 2.5%

Processing Features

Patient samples (100 µL/well): ~16 minutes/full plate
 Reagent (100 µL/well): ~4 minutes/full plate

Data Processing

User Interface
 Software environment: Windows XP operating system

Communication

RS-232 Serial Port
 On-line (bi-directional): ASTM format

Dimensions

Processor module
 Size in inches (H x W x D): 38.0" x 45.2" x 30.4"
 Weight: 209 lbs.

Table Module

Size in inches (H x W x D): 22.0" x 48.0" x 32.0"

Electrical

Line voltage (adjustable): 110/260 V
 Frequency: 47-63 Hz
 Power consumption: 500 VA

Internal Bar Code Scanner

Positive sample ID: yes
 Reagent ID: yes
 Reads: Interleaved 2 of 5, Code 3 of 9, IATA 2 of 5, Industrial 2 of 5, UPC A & E, EAN 8 or 13, Code 128, Pharmacode, Codabar & EAN Addendum 2 or 5

Note: Specifications are subject to change without prior notice. EVOLIS is a trademark of Bio-Rad Laboratories, Inc. Windows is a trademark of Microsoft Corporation.

Ordering Information

Catalog No.	Description	
89700J	EVOLIS™ System (including analyzer, PC, monitor and printer).....	1 system
25167	EVOLIS™ Workstation Table	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

Clinical
 Diagnostics Group

Website 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