# FAST... LOW SAMPLE VOLUME... RELIABLE VI-CELL MetaFLEX ANALYZER









# **Fast**

Results in 35 sec.

# **Small Sample Volume**

65µL sample volume

# Reliable

Ready for use when you need it, high uptime of 23 hours per day

# R&D, QC and Manufacturing Applications

Ideal for micro to large scale cell culture applications, the Vi-CELL MetaFLEX is designed for rapid and accurate Bioanalyte analysis.

# Analyzer Performance

Measuring System	
Sample volume (all parameters)	65 µL
Measuring time (all parameters) 35 sec*	
Cycle time	60 sec*
Throughput	44 samples / per hour*
Average uptime	23.5 hours / per day**

<sup>\*</sup> May vary during startup

<sup>&</sup>quot;May vary during startup; 2.5 min for system calibration is the longest activity

Sensor Cassette	
In-use lifetime	30 days (or until max number of tests)
Shelf life	4 months
Storage temperature	2-8° C / 35-46° F
Automatic QC	Yes
Number of tests	300

Solution Pack	
In-use lifetime	30 days (or until solution pack is empty)
Shelf life	6 months
Storage temperature	2-25° C / 35-77° F

# **Quality Management**

## Automatic quality management system

- 3 dedicated QC solutions
- · Automatic detection and correction of failures
- Continuous system and analysis checks
- Air detection
- · Automatic lockout of parameter that fails QC
- Customizable QC schedule

Measured Parameters				
Туре	Parameter	Units	Range of indication	
рН	На	pH scale	6.3 - 8.0	
Cell Culture Gas	pCO <sub>2</sub>	mmHg; Torr	5 - 250	
		kPa	0.67 - 33.3	
		mmHg; Torr	0 - 803	
	pO <sub>2</sub>	kPa	0 - 107	
Electrolyte	cK+	mmol/L	0.5 - 25.0	
		meq/L	0.5 - 25.0	
	cNa⁺	mmol/L	7 - 350	
	CING	meq/L	7 - 350	
		mmol/L	<0.1 - 9.99	
	cCa²⁺	meq/L	<0.1 - 19.98	
		mg/dL	<0.1 - 40.04	
	cCl <sup>-</sup>	mmol/L	<0.1 - 350	
		meq/L	<0.1 - 350	
Metabolite		mmol/L	0 - 60	
	<i>c</i> Glu	mg/dL	0 - 1081	
		g/L	0 - 10.81	
		mmol/L	O - 31	
	cLac	meq/L	O - 31	
		mg/dL	0 - 279	
		g/L	0 - 2.79	

The Range of indication for a parameter is the range within which the analyzer is physically capable of measuring.

# Ordering Information

Part Number	Description
B93015	Vi-CELL MetaFLEX
B93029	Vi-CELL MetaFLEX Solution Pack
B93030	Vi-CELL MetaFLEX Sensor Cassette



# Hardware

### Computer specifications

- Processor Intel Celeron® M 600 MHz with 512K L2 Cache
- 1 GB RAM
- 2 GB SolidState storage
- 8.4" color TFT-LCD, resolution 800 × 600 Touch screen
- 4" thermal-sensitive printer

- Left/right hand operation
- Position for syringe as well as capillary and test tube
- Aspiration from capillary without adapter
- Aspiration time 5 seconds

### Interface

- Built-in barcode reader for operator & sampler ID Accepted codes: UPC/EAN, Code 128, Code 39, Code 93, I 2 of 5, Discrete 2 of 5, Codabar and more
- Serial interface RS232 with power for external barcode reader
- 3 USB connections
- · Optional external keyboard
- Optional external mouse
- Optional external barcode reader

# Software

### Software platform

- Windows® 10 Enterprise IoT
- Sybase® SQL Anywhere

### Data capacity

- · Sample results log: 2000
- · Activity log: 5000
- · Calibration adjustment log: 1000
- · Data secured by password protection
- · 8 different user profiles
- Unlimited ID access verification
- · Printer display options
- Auto print (on/off)
- Select derived parameters
- Select input variables
- Reference ranges with results

### Security and QA features

- Advanced planning of replacement and QC schedules
- Optional automatic QC at startup and after replacements
- Customizable QC and calibration schedule

### Communication

- · HIS/LIS communication
- High-level protocols:
  - ASTM
  - HI 7
  - POCT1-A
- Low-level serial protocols:
  - ASTM 1381-91
  - E1394-91
  - Serial RAW
- · Low-level network protocols:
  - TCP/IP

# Additional Information

### Dimensions

45 cm 17.7 in Height 25 cm 9.8 in Width Depth 29 cm 11.4 in Weight 11 kg 24.4 lbs

### Other

· Startup time Operating

environment

correction

Power Thermostat control

Altitude

Up to 4000 m / 13,124 feet above sea level

1.5 hours average (less than 4 hours)

15 - 32° C / 59 - 89.6° F, 20 - 80% RH

100 - 240 V; 50 - 60 Hz; 250 VAC Sensor cassette: 37 ± 0.15° C /

98.6 ± 32.27° F Oximetry: 37 ± 0.30° C / 98.6 ± 32.54° F

CLASS 1 LASER PRODUCT

The Vi-CELL MetaFLEX ANALYZER and reagents are not intended or validated for use in the diagnosis of disease or other conditions.

© 2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks used herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.



For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com