



# Fast, Reliable & Low-Volume Analysis of Bioanalytes

Vi-CELL MetaFLEX Analyzer

ACCELERATING  
*answers*

 **BECKMAN  
COULTER**  
*Life Sciences*

# Rapid & Accurate Bioanalyte Analysis for R&D, QC & Manufacturing Applications



## Vi-CELL MetaFLEX Analyzer performance

### Measuring System

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

\* May vary during startup

\*\* 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	60 days
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	60 days
Storage temperature	2-25° C / 35-77° F



### Fast

Results in 35 sec.



### Small sample volume

65 µL sample volume



### Reliable

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

## Automatic quality management system provides:

- 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

Type	Parameter	Units	Range of indication
pH	pH	pH scale	6.3 - 8.0
Cell Culture Gas	pCO <sub>2</sub>	mmHg; Torr	5 - 250
		kPa	0.67 - 33.3
	pO <sub>2</sub>	mmHg; Torr	0 - 803
		kPa	0 - 107
Electrolyte	cK <sup>+</sup>	mmol/L	0.5 - 25.0
		meq/L	0.5 - 25.0
	cNa <sup>+</sup>	mmol/L	7 - 350
		meq/L	7 - 350
	cCa <sup>2+</sup>	mmol/L	<0.1 - 9.99
		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	cGlu	mmol/L	0 - 60
		mg/dL	0 - 1081
		g/L	0 - 10.81
	cLac	mmol/L	0 - 31
		meq/L	0 - 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 Analyzer
B93029	Vi-CELL MetaFLEX Solution Pack
B93030	Vi-CELL MetaFLEX Sensor Cassette



# Vi-CELL MetaFLEX Specifications

## Hardware

### Computer specifications

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

### Inlet

- 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, 1 2 of 5, Discrete 2 of 5, Codabar and more
- Serial interface RS232 with power for external barcode reader
- 3 USB connections
- Optional external devices:
  - Keyboard
  - Mouse
  - Barcode reader

## Software

### Software platform

- Windows® 10 Enterprise IoT
- Sybase® SQL Anywhere

### Data capacity

- Sample results log: 2,000
- Activity log: 5,000
- Calibration adjustment log: 1,000
- 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
  - HL7
  - POCT1-A
- Low-level serial protocols:
  - ASTM 1381-91
  - E1394-91
  - Serial RAW
- Low-level network protocols:
  - TCP/IP

## Additional Information

### Dimensions

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

### Other

- Startup time 1.5 hours average (< 4 hours)
- Operating environment 15 - 32° C / 59 - 89.6° F, 20 - 80% RH
- Altitude correction Up to 4,000 m / 13,124 feet above sea level
- Power 100 - 240 V; 50 - 60 Hz; 250 VAC
- Thermostat control Sensor cassette: 37 ± 0.15° C / 98.6 ± 32.27° F  
Oximetry: 37 ± 0.30° C / 98.6 ± 32.54° F



**Learn more**  
about the Vi-CELL  
MetaFLEX Analyzer



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.

© 2026 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)  
2023-GBL-EN-103709-v2

