

## It's not just a UV/Vis. It's a LAMBDA.

### The Bio-calculator – with the heart of a LAMBDA

Introducing the PerkinElmer LAMBDA™ Bio- and Bio+ – personal UV/Vis spectrophotometers for all your routine life-science calculations. Continuing our tradition of providing you with quality UV/Vis instruments for more than 60 years, these low-cost spectrophotometers have been designed to acquire and calculate results in a matter of seconds. With the same simplicity as a desktop calculator, the LAMBDA Bio and Bio+ have a series of preprogrammed methods to support the demands of both molecular biology and biotech laboratories:

- DNA, RNA, concentration and purity, and Oligonucleotide methods
- Protein methods, including: Direct UV, BCA, Bradford, Lowry, Buiret, and Cell-density measurements
- Curve fitting with up to nine standards in triplicate assures accurate results

### Performance that wow's the crowd and the finance people

With its large LCD screen and full scanning capabilities, the LAMBDA Bio is a crowd pleaser. Built to last, easy to use, and even easier to own. No moving parts, virtually no maintenance, and no consumables make the LAMBDA Bio the perfect tool when the instrument just needs to work, day-in and day-out. Rely on the LAMBDA Bio. It's the smart choice for your laboratory in any application, from teaching – where operator training may be an issue, to the most advanced R&D laboratory – where precious samples need to be processed right, the first time. The LAMBDA Bio will win the hearts and minds of anyone working with it. After all, it is a LAMBDA.



LAMBDA Bio and Bio+ Spectrophotometer



## **Designed To Make Spectrophotometry Easier**

# You don't need an impressive budget to have an impressive instrument

The performance you'll experience with the LAMBDA Bio and Bio+ is only normally associated with much more expensive designs. Experience:

- Instant answers Unlike other instruments on the market, our LAMBDA Bio requires no warm-up time. Our ultra-fast initialization will allow you to capture a full range of spectrum in less than three seconds.
- **Continuous uptime** Our maintenance-free design with no moving parts and Xenon light source provides you with the reliability your laboratory needs.
- **Superb stability** Avoid instrument drift from excessive heat with our split-beam design.
- **Ruggedness** Our sealed, zero-bounce membrane keypad is resistant to even the largest solvent spills.

#### Provides you with the versatility to do just about anything

Your work environment may be complex but that doesn't mean your instrument needs to be. With the simplicity and elegance of the LAMBDA Bio, you'll be empowered to analyze samples when and where your projects demand:

- View, expand, and manipulate full spectra quickly and easily with the clear 11.5 cm by 8.7 cm backlit LCD display.
- Feel confident Low-stray light provides exceptional linearity and confidence in your measurements.
- **Keep up the pace** Our innovative polychromator technology will give you the speed and performance you need to keep up with your fast-paced laboratory.
- **Save time** Our optical design allows samples to be run with no sample compartment lid.
- **Personalize your laboratory** The huge on-board memory will allow you to store up to 90 methods along with unique method names and calibration information (only available on the Bio+).
- **Convenient sampling flexibility** Measure your samples in any of our high-quality, precision glass, quartz, or disposable cells:
  - Macro cells: 3.5 ml
     Semi-micro cells: 1.4 ml
     Ultra-micro cells: 15 μl



## Share the knowledge – multiple archiving and data sharing options

- Collect data from up to eight instruments A single PC can be configured to accept and automatically display data from up to eight instruments. Data can be archived in Excel® (.xls), ASCII (.csv or .txt), graphics (.emf), or rich text (.rtf) formats.
- Built-in USB port Transfer data to your computer for viewing, manipulating, printing, or archiving using the built-in USB connection port. LAMBDA Bio PVC software is provided as standard with all instruments.
- Optional integrated thermal printer For your permanent records, use the optional thermal printer to print all of your results. With the standard 57 mm printer roll, your print-out can accommodate Calibration Data, Result, and Time and Date.
- Secure Digital (SD) Flash Storage Card or Bluetooth®
   compatible For your archiving convenience, the LAMBDA Bio
   can be outfitted to accommodate a standard SD flash storage
   card or Bluetooth® accessory allowing you additional flexibility
   for transferring data.

#### Experience the power - flexible on-board software

The LAMBDA Bio software is optimized to provide you with answers fast and conveniently. Save your own methods or choose from our broad range of pre-loaded measurement modes for simple, efficient analysis.

Calculates cell density to ensure cell proliferation is sufficient for induction or harvesting. Choose from OD units or enter a factor to calculate cells/ml.



Sample	0.189 A	A230
6	0.390 A	A260
	0.009 A	A280
Concentration	0.000 A	A320
19.50	A280	A260
	43.3	
Units	A260/A230	
µg/ml	2.063	

Measure DNA, RNA, or Oliglo-
nucleotides. Pathlength and
dilution factors produce direct
concentration results.

A260	Sample
0.495 A	7
A280	Result
0.036 A	000
A320	-322
0.002 A	
	Units
	ug/ml

Calculate protein concentration by direct UV, BCA, Biuret, Lowry, and Bradford methods.

	S	ample
		1
	λ	620 nm
		Result
$   $ $\wedge$ $ $	U.	038
1/ 1/		

Allows you to specify a single wavelength to acquire data in %T or absorbance.

Wavelength	Sample
261 nm	7
Absorbance	Concentration
0.016 A	0.800
Factor	0.000
50.0	
	Units
	μg/ml

Data acquired at a defined wavelength will be multiplied by a constant selected by the user and the result displayed.

Specifications		LAMBDA Bio	LAMBDA Bio+
Optical design		Split beam	
Wavelength	Range Scanning range Calibration Accuracy Reproducibility	190-1100 nm 200-950 nm Automatic when the instrument is switched on ±2 nm ±1 nm	
Spectral bandwidt	h	5 nm	3 nm
Light source		Pulsed Xenon lamp – 5+ year lifetime	
Photometric Range Linearity  Reproducibility		-0.300 to 2.500 A, 0 to 199 %T ±0.005 A or 1% of the reading, whichever is the greater at 546 nm ±0.003 A (0-0.5 A), ±0.007 A (0.5-1.0 A)	
Stray light		<0.5% at 220 nm and 340 nm using NaNO <sub>2</sub>	
Zero stability		±0.01 A/hour	
Noise		0.005 peak to peak 0.002 rms	
Stored capacity		9	90
Display		11.5 cm x 8.8 cm	
Keypad		Sealed membrane keypad	
Outputs		USB, SD Card (optional), Bluetooth® (optional), Printer (optional)	
Dimensions (WxHx	(D)	340 mm x 330 mm x 170 mm (340 mm x 420 mm x 170 mm with printer attached)	
Weight		3.2 kg (7.1 lb)	
Power input		100-250 V, 50/60 Hz, Max 30 VA	
Languages		English, Spanish, German, French, Italian, Japanese	
Warranty		1 year	

Online Ordering Information		Options and Consumables		
Part No.	System	Part No.	Description	
L7110184	LAMBDA Bio	L7110230	Printer module	
L7110185	LAMBDA Bio with Printer	L7110231	PVC Software and PC USB Connection*	
L7110186	LAMBDA Bio+	L7110232	Spare printer paper (20 rolls)	
L7110187	LAMBDA Bio+ with Printer	L7110233	Bluetooth® accessory	
L7110188	LAMBDA Bio+ with SD Card	* Included in	all models except the LAMBDA Bio and LAMBDA Bio with printer.	

### Order online at www.perkinelmer.com/lambdabio

PerkinElmer, Inc. 940 Winter Street Waltham, MA 02451 USA P: (800) 762-4000 or (+1) 203-925-4602 www.perkinelmer.com

