

Personal UV/Vis Spectrophotometers



New Personal Life-Science UV/Vis Spectrophotometers – Speed, Stability and Confidence Results

- No moving parts for the ultimate in stability and reliability
- Xenon source – typical lifetime of five years
- Compact design uses minimal bench space
- Scans full spectra almost instantaneously
- Transfer data to your computer using an SD memory card or USB PC connector
- Large, intuitive full graphics display
- Optional integrated thermal printer
- Multiple language operation as standard

LAMBDA BIO AND BIO+

The LAMBDA™ Bio and Bio+ have been specially designed for life-science laboratories that require a dedicated instrument for the determination of nucleic acid purity and concentration, protein concentrations and cell density measurements. The LAMBDA Bio and Bio+ are compact, lightweight and designed for maximum convenience and ease-of-use.

- Full suite of stored life-science methods
- Compatible with low volume cuvettes

LAMBDA BIO

a personal spectrophotometer for nucleic acid and protein determinations, cell density measurements and more

LAMBDA BIO+

all the features of the LAMBDA Bio plus routing assays, kinetics, rapid scanning and concentration measurements

Specifications	LAMBDA Bio and Bio+
Wavelength Range	190 - 1100 nm
Absorbance Range	-0.3 to 2.5 A
Bandwidth	5 nm (3 nm LAMBDA Bio+)
Stray Light	0.5% T at 220 nm and 340 nm
Wavelength Accuracy	±2 nm
Photometric Reproducibility	±0.002A at 0 - 0.5A, 546 nm
Stored Methods	9 (90 LAMBDA Bio+)
Outputs	USB, Printer (option), Additional SD Card (option), Bluetooth (option) on LAMBDA Bio+
Dimensions (W x D x H)	340 mm x 330 mm x 170 mm
Weight	<3.2 kg

Ordering Information

Part No.	System
L7110184	LAMBDA Bio
L7110185	LAMBDA Bio with Printer
L7110186	LAMBDA Bio+
L7110187	LAMBDA Bio+ with Printer
L7110188	LAMBDA Bio+ with SD Card

Options and Consumables

Part No.	Description
L7110230	Printer module
L7110231	Print utility and cable**
L7110232	Spare printer paper (20 rolls)
L7110233	Bluetooth accessory

** Included as standard with the LAMBDA Bio+

LAMBDA XLS AND LAMBDA XLS+

The LAMBDA XLS and XLS+ have been specially designed for academic and industrial laboratories that require a small-footprint, dedicated instrument for a wide range of applications. The LAMBDA XLS and XLS+ are extremely reliable and sensitive instruments that provide a level of performance which is only normally found in more expensive systems.

- Scans full spectra almost instantaneously
- Measure using 10 mm, 20 mm or 40 mm pathlength cells

Specifications	LAMBDA XLS and XLS+
Wavelength Range	190 - 1100 nm
Absorbance Range	-0.3 to 2.5 A
Bandwidth	5 nm (3 nm LAMBDA XLS+)
Stray Light	0.5% T at 220 nm and 340 nm
Wavelength Accuracy	±2 nm
Photometric Reproducibility	±0.002A at 0 - 0.5A, 546 nm
Stored Methods	90
Outputs	USB, SD Card (option), Bluetooth (option), Printer (option)
Dimensions (W x D x H)	340 mm x 330 mm x 170 mm
Weight	<3.2 kg

LAMBDA XLS

a personal 5 nm spectrophotometer to meet your scanning and fixed wavelength studies, kinetics and concentration measurements needs

LAMBDA XLS+

a higher resolution (3 nm bandwidth) version of the LAMBDA XLS

Ordering Information

Part No.	System
L7110189	LAMBDA XLS
L7110190	LAMBDA XLS with Printer
L7110191	LAMBDA XLS with SD Card
L7110192	LAMBDA XLS+
L7110193	LAMBDA XLS+ with Printer
L7110194	LAMBDA XLS+ with SD Card

Options and Consumables

Part No.	Description
L7110230	Printer module
L7110231	Print utility and cable
L7110232	Spare printer paper (20 rolls)
L7110233	Bluetooth accessory