

UV/VIS  
UV/VIS/NIR Spectroscopy**Features and Benefits**

- Large, flexible sampling capacity
- Fast, accurate random access
- Corrosion-resistant sampling components
- Automatic rinsing
- Flow through rinse station to minimize sample-to-sample contamination

## S10 Autosampler for UV/VIS, UV/VIS/NIR Spectroscopy

The PerkinElmer S10 autosampler is a computer-controlled, multi-purpose sampling system for the Lambda UV/VIS, UV/VIS/NIR spectrophotometers.

The autosampler automates standard and sample introduction for instrument

calibration and sample analysis, extending the spectrometer's capabilities to those of a fully automated analytical workstation.

**Fast, accurate sample access**

An advanced drive system with high-precision motors moves the sampling arm of the S10 in the X and Y coordinates simultaneously, minimizing changeover time between samples. Random-access programming gives the user exceptional flexibility in the placement of samples and reference solutions.

**Large, flexible sample capacity**

The S10 provides more than 200 vessel positions for maximum capacity. Sample racks are easily interchanged, reducing analysis setup times and giving you the flexibility to use different-sized sample vessels matched to your analytical needs.

The S10 autosampler has a built-in peristaltic pump for continuous rinsing of the sampling capillary between samples to significantly reduce the risk of carryover. The sample probe, including the sample transport capillary tubing, is acid- and solvent-resistant for long-term operation.

## Specifications

### S10 Autosampler for UV/VIS Part No. N2020004

#### Sampling Arm Movement

Maximum movement	Vertical: 145 mm; Horizontal: x 302 mm, y 222 mm
Resolution	0.1 mm $\pm$ 4%
Speed (movement from one sampling position to the next)	Typically 3 seconds or less

#### Power Requirements

S10	Built-in power supply: 115/230 V, 50/60 Hz.
-----	---

#### Dimensions

Width x Height x Depth	46 cm x 53 cm x 38 cm
Weight	5 kg

## Rack Configurations

#### Tray Type

#### Required Sub Trays

Tray E	1 tray of B3140621 and 2 trays of B3140617
Tray F	1 tray of B3001647 and 2 trays of B3140618
Tray G	3 trays of B3140621

## Sample Trays

Order No.	B3140617	B3140618	B3140621	B3001647
Sample locations	90	60	21	9/29
Location diameter (mm)	13	16	30	30/16
Sample vessel (mL)	6 or 8	15	50	50/15

## Polypropylene Tubes

We offer a wide variety of high-quality, capped and uncapped autosampler tubes and supplies for your S10 Autosampler. Ranging in size from 8 mL to 50 mL, these tubes are available with either round or conical bottoms. Free-standing options are also available.\*



#### Autosampler Tubes

Part No.	Capacity	Size	Type	Qty
N0777156	8 mL	13 mm x 100 mm	Round Bottom	1200
N0777167	15 mL	17 mm x 100 mm	Round Bottom	1200
N0777599		Plug for 15 mL Vials (N0777167)		1200
N9301205	16 mL	17 mm x 100 mm	Round Bottom	1000
N0777158	50 mL	30 mm x 115 mm	Round Bottom	500

#### Autosampler Tubes with Caps

Part No.	Capacity	Size	Type	Qty
B0193233	15 mL	13 mm x 100 mm	Conical Bottom	500
B0193234*	50 mL	30 mm x 115 mm	Conical Bottom	500

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2007-2015, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.