

Atomic Spectroscopy

Key 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 Atomic Spectroscopy

The PerkinElmer® S10 autosampler is a computer-controlled, multi-purpose sampling system for AA, ICP-OES and ICP-MS. 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

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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2007-2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

007899B_01