

PINAACLE D900 AA CONSUMABLES AND SUPPLIES



INTRODUCTION AND OVERVIEW

PinAAcle™ D900 Atomic Absorption (AA) consumables and supplies are designed with your instrument in mind. Each part is fit to perform and manufactured with the highest quality materials available. Whether you are looking for lamps, graphite tubes, sample introduction components or standards, we have the consumables you need to keep your PinAAcle D900 AA spectrometer up and running smoothly and efficiently.

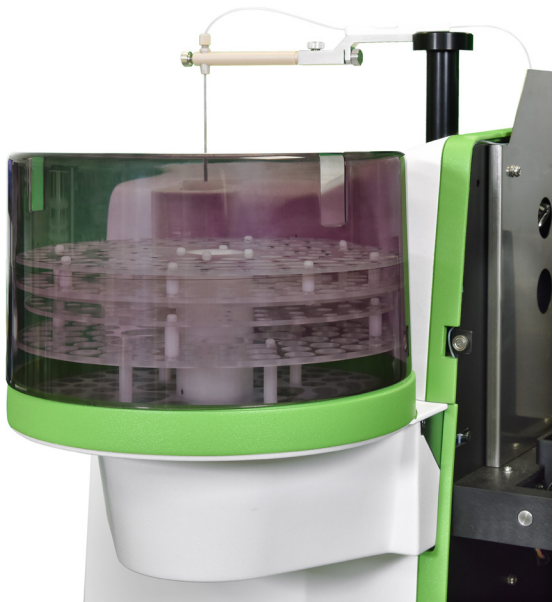
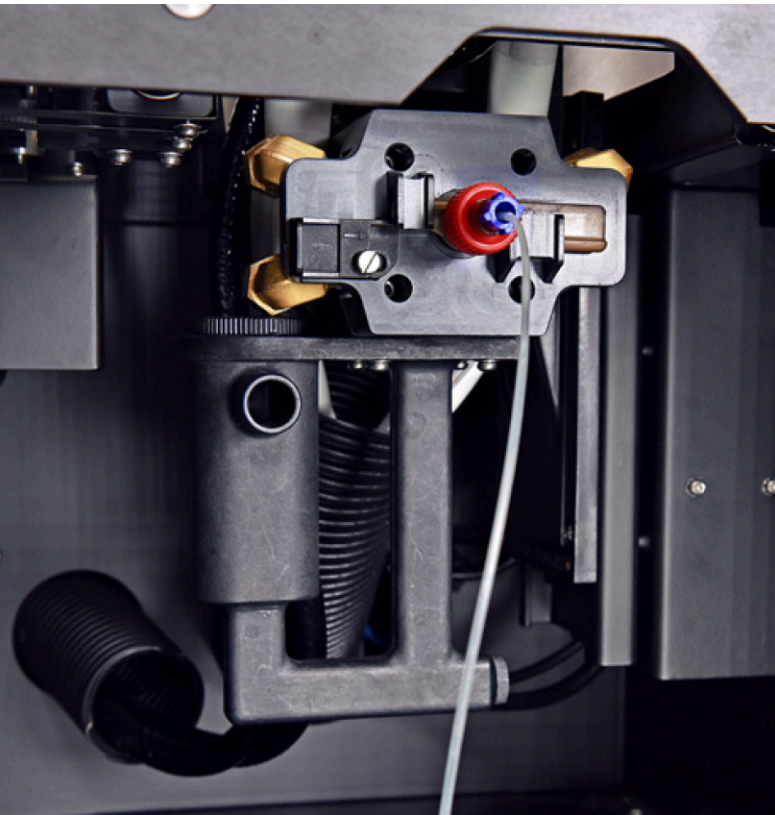


TABLE OF CONTENTS



Hollow Cathode Lamps.....	3
Deuterium Lamps	4
HGA Graphite Tubes.....	4
Contact Cylinders	4
Graphite Furnace Autosampler Cups	4
Autosampler Tubes	4
Burner Heads.....	4
Nebulizers.....	4
Sampler Probes and Spares	5
Heating Wire	5
N ₂ O Hose Assembly.....	5
Matrix Modifiers.....	5

TRUST THE CONSUMABLES
ENGINEERED FOR YOUR INSTRUMENTS

TRUST THE CONSUMABLES ENGINEERED FOR YOUR INSTRUMENTS

Single-Element Hollow Cathode Lamps

PerkinElmer is the world leader in atomic absorption. Atomax™ 1.5 inch coded hollow cathode lamps are crafted to the same high standards that have made our genuine PerkinElmer 2 inch Lumina™ lamps the best in the industry. Drawing on more than 40 years of HCL design and manufacturing expertise, you can trust the consumables that were engineered for your instrument.

Single-Element Atomax Coded Hollow Cathode Lamps		
Element Name	Symbol	Part No.
Aluminum	Al	N3300603
Antimony	Sb	N3300661
Arsenic	As	N3300615
Barium	Ba	N3300616
Beryllium	Be	N3300619
Bismuth	Bi	N3300628
Cadmium	Cd	N3300601
Calcium	Ca	N3300610
Cesium	Cs	N3300617
Chromium	Cr	N3300606
Cobalt	Co	N3300620
Copper	Cu	N3300604
Gold	Au	N3300662
Iron	Fe	N3300612
Lead	Pb	N3300602
Lithium	Li	N3300630
Palladium	Pd	N3300624
Platinum	Pt	N3300664
Magnesium	Mg	N3300611
Manganese	Mn	N3300613
Molybdenum	Mo	N3300627
Mercury	Hg	N3300614
Nickel	Ni	N3300605
Phosphorus	P	N3300667
Potassium	K	N3300609



Atomax Coded Lamp

Single-Element Atomax Coded Hollow Cathode Lamps		
Element Name	Symbol	Part No.
Selenium	Se	N3300666
Silicon	Si	N3300665
Silver	Ag	N3300623
Sodium	Na	N3300608
Tin	Sn	N3300629
Titanium	Ti	N3300621
Thallium	Tl	N3300625
Vanadium	V	N3300618
Zinc	Zn	N3300607
Antimony	Sb	N3300670
Arsenic	As HPHCL	N3300626
Cadmium	Cd HPHCL	N3300622
Selenium	Se HPHCL	N3300669
Tin	Sn HPHCL	N3300668

Background Corrector Source	
Description	Part No.
Deuterium Lamp	09201018



Deuterium Lamp

HGA Graphite Tubes

Description	Part No.
HGA Graphite Tubes with Integrated Platform Pkg. 3	N3303510
HGA Graphite Tubes with Integrated Platform Pkg. 5	N3303508
HGA Graphite Tubes with Integrated Platform Pkg. 10	N3303509



HGA Graphite Tubes

Contact Cylinders

Description	Part No.
Contact Cylinders 1-Set	B0128495
Graphite Guide Tool	N3302530



Contact Cylinders

Graphite Furnace Autosampler Cups

Volume	Composition	Qty.	Part No.
2.5 mL	Polypropylene	1,000	B3001566
7.0 mL	Polypropylene	100	B3001567



Graphite Furnace Cups

Autosampler Tubes

Description	Size	Qty.	Part No.
15 mL Round Bottom Tubes without Plugs	17 mm x 100 mm	1,200	N0777167
15 mL Conical Tubes with Red Flat Caps	17 mm x 120 mm	500	B0193233
50 mL Freestanding Tubes with Red Flat Caps Attached	28 mm x 115 mm	500	B0193234



Autosampler Tubes

Burner Heads

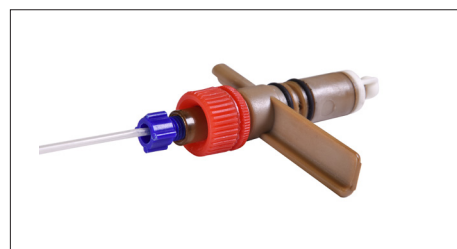
Description	Part No.
Burner Head Main Body	N3306712
Burner Head N ₂ O Assembly	N3306718
Air Burner Head	N3308731
N ₂ O ID Tip Burner Head	N3308733
Burner Head O-Ring 15.54 mm x 2.62 mm	09220611



Burner Head

Nebulizers

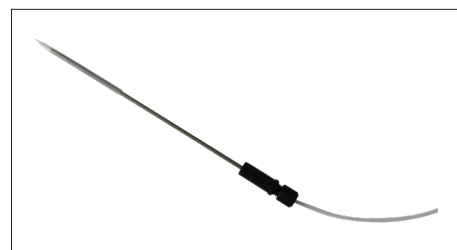
Description	Part No.
Nebulizer Assembly	N3160144



Nebulizer Assembly

Sampler Probes and Spares

Description	Part No.
Sampler Probe Furnace	N3306721
Sampler Probe Flame	N3306722
O-Ring 1.42 mm ID x 1.52	09200960
Autosampler Rinse Outer Tube	N3302833
Furnace Autosampler Sample Capillary	N3302837
Flame Autosampler Sample Capillary	N3302838
Peristaltic Pump Inner Tube	N3302841



Sampler Probe

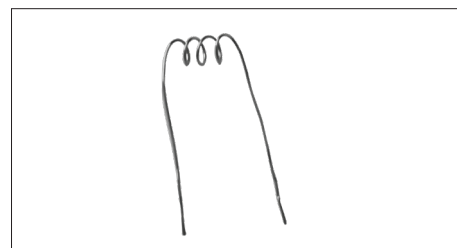
N₂O Hose Assembly

Description	Part No.
N ₂ O Hose Assembly	N3150176

N₂O Hose Assembly

Heating Wire

Description	Part No.
Heating Wire	N3303550



Heating Wire

Matrix Modifiers

Modifiers	Composition	Qty.	Part No.
Mg(NO ₃) ₂	1% Mg (as Nitrate)	100 mL	B0190634
Pd	1% Pd (as Nitrate)	50 mL	B0190635
NH ₄ H ₂ PO ₄	10% NH ₄ H ₂ PO ₄	50 mL	N9303445



Matrix Modifiers

Everyday you count on PerkinElmer to provide you with solutions that deliver reliable performance, control operating costs and maximize operational time. Our complete portfolio of consumables, parts, supplies, training and service helps you meet both routine and demanding measurement challenges. We invest heavily in testing and validating our products to ensure you receive guaranteed compatibility and performance – on-time, for every instrument in your laboratory.

Shop Online Today!

www.perkinelmer.com/buy

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 ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.