



AA Graphite and Lamps for Non-PerkinElmer Instruments

Atomic Absorption Graphite Furnace Tubes

PerkinElmer supplies a wide range of graphite tubes for non-PerkinElmer instrumentation. Graphite tubes are made with the highest quality material guaranteed to work for your AA instrument.

Use the convenient tables below to identify the compatible graphite furnace tube for your AA instrument.

GBC Instruments

Part No.	GBC Part No.	Description	Qty./pack
N9304600	99-0059-00	Standard tube, pyrolytically coated	10
N9304601	99-0060-00	Platform, pyrolytically coated	10
N9304602	99-0342-00	Standard tube, coated, with Omega-platform	10
N9304603	99-0061-00	Electrode contacts	2
N9304604	45-0004-00	Shroud	1

Hitachi® Instruments

Part No.	Hitachi® Part No.	Description	Qty./pack
N9304605	180-7444	Standard tube, pyrolytically coated	10
N9304606	180-7400	Standard tube, uncoated	10
N9304607	180-7404	Standard platform, pyrolytically coated	10
N9304608	190-6003	Tube for extended injection volume, pyrolytically coated	10
N9304609	180-7401	Contacts, AD 20 mm	2
N9304610	172-8805	Zeeman contacts, AD 18 mm	2
N9304611	-	Omega Tube	10

Shimadzu® Instruments

Part No.	Shimadzu® Part No.	Description	Qty./pack
N9304612	206-50588	Standard tube, pyrolytically coated (90° contact cone)	10
N9304613	206-50587	Standard tube, uncoated (90° contact cone)	10
N9304614	206-5887-02RI	Tube with built-in platform, coated (90° contact cone)	10
N9304615	206-50602	Left Electrode with housing (90° cone)	1
N9304616	206-50603	Right Electrode (90° cone)	1

Thermo™ Instruments

Part No.	Thermo™ Part No.	Description	Qty./pack
N9304617	9423 393 95031	Partridged tube, uncoated	10
N9304618	9423 393 95071	Partridged tube, coated	10
N9304619	9423 393 95041	Partridged tube, coated (extended lifetime)	10
N9304620	9423 393 95161	Zeeman Electrode Contact	2
N9304621	9423 393 95011	Standard Electrode Contact	2
N9304622	9423 490 20101	Omega Tube	10

Varian™ Instruments			
Part No.	Varian™ Part No.	Description	Qty./pack
N9304623	63-100011-00	Plateau tube, pyrolytically coated	10
N9304624	63-100013-00	Bone platform, pyrolytically coated (for use in plateau tubes)	10
N9304625	63-100012-00	Partition tube, pyrolytically coated	10
N9304626	63-100015-00	Partition tube, uncoated	10
N9304627	63-100012-EL	Partition tube, pyrolytically coated extended life	10
N9304628	63-100037-00	Omega-platform tube	10
N9304629	63-100018-00	Shroud	1
N9304630	63-100016-00	Electrode	2
N9304631	63-100019-00	Zeeman Shroud	1
N9304632	63-100017-00	Zeeman Electrode (Older Instruments)	2
N9304633	63-100035-00	Zeeman Electrode (Newer Instruments)	2
N9304634	63-100036-00	Zeeman Shroud GTA120, pyrocoated	1

Atomic Absorption 1.5" Hollow Cathode Lamps

PerkinElmer supplies a wide range of single-element hollow cathode lamps for non-PerkinElmer instrumentation. HCL lamps are made with the highest quality of materials guaranteed to work for your AA instrument. Atomax™ 1.5" hollow cathode lamps are crafted to the same high standards that have made PerkinElmer 2" Lumina™ hollow cathode lamps the best in the industry. Drawing on more than 40 years of HCL design and manufacturing expertise, Atomax lamps now offer laboratories around the world a dependable, top-quality lighting source, regardless of the brand of AA instruments used.

Use the convenient table below to identify the compatible Atomax lamp for your AA instrument.

Element Name	Symbol	PerkinElmer Part No.	Varian™ Part No.	Cathodeon Part No.	Shimadzu® Part No.	Thermo™ Part No.	Photron Part No.
Aluminum	Al	N2025301	5610122000	3UNX/Al	200-38422-01	9423 390 20131	P801
Antimony	Sb	N2025347	5610122100	3QNY/Sb	200-38422-24	9423 390 20511	P802
Arsenic	As	N2025302	5610122200	3QNY/As	200-38422-42	9423 390 20331	P803
Barium	Ba	N2025305	5610122300	3BAX/Ba	200-38422-03	9423 390 20561	P804
Beryllium	Be	N2025306	5610122400	3QNY/Be	200-38422-04	9423 390 20041	P805
Bismuth	Bi	N2025307	5610122500	3QNY/Bi	200-38422-43	9423 390 20831	P806
Boron	B	N2025304	5610122600	3UNX/B	200-38422-39	9423 390 20051	P807
Cadmium	Cd	N2025309	5610122700	3QNY/Cd	200-38422-06	9423 390 20481	P808
Calcium	Ca	N2025308	5610122900	3UNX/Ca	200-38422-05	9423 390 20201	P809
Cesium	Cs	N2025312	5610122800	3BNX/Cs	200-38422-27	9423 390 20551	P810
Chromium	Cr	N2025311	5610123100	3UNX/Cr	200-38422-07	9423 390 20241	P812
Cobalt	Co	N2025310	5610123200	3UNX/Co	200-38422-09	9423 390 20271	P813
Copper	Cu	N2025313	5610123300	3UNX/Cu	200-38422-08	9423 390 20291	P814
Dysprosium	Dy	N2025314	5610123400	3BNX/Dy	200-38422-60	9423 390 20661	P815
Erbium	Er	N2025315	5610123500	3BNX/Er	200-38422-61	9423 390 20681	P816
Europium	Eu	N2025316	5610123600	3BAX/Eu	200-38422-62	9423 390 20631	P817
Gadolinium	Gd	N2025319	5610123700	3UNX/Gd	200-38422-63	9423 390 20641	P818
Gallium	Ga	N2025318	5610123800	3UNX/Ga	200-38422-40	9423 390 20311	P819
Germanium	Ge	N2025320	5610123900	3UNX/Ge	200-38422-11	9423 390 20321	P820
Gold	Au	N2025303	5610124000	3UNX/Au	200-38422-25	9423 390 20791	P821
Hafnium	Hf	N2025321	5610124100	3UAX/Hf	200-38422-64	9423 390 20721	P822
Holmium	Ho	N2025323	5610124200	3BNX/Ho	200-38422-65	9423 390 20671	P823
Indium	In	N2025324	5610124400	3UNX/In	200-38422-48	9423 390 20491	P824
Iridium	Ir	N2025325	5610124500	3QNY/Ir	200-38422-48	9423 390 20771	P825

PerkinElmer Atomax

Element Name	Symbol	PerkinElmer	Varian™	Cathodeon	Shimadzu®	Thermo™	Photron
		Part No.	Part No.	Part No.	Part No.	Part No.	Part No.
Iron	Fe	N2025317	5610124600	3UNX/Fe	200-38422-10	9423 390 20261	P826
Lanthanum	La	N2025327	5610124700	3UAX/La	200-38422-29	9423 390 20571	P827
Lead	Pb	N2025339	5610124800	3QNY/Pb	200-38422-21	9423 390 20821	P828
Lithium	Li	N2025328	5610124900	3BAX/Li	200-38422-30	9423 390 20031	P829
Lutetium	Lu	N2025329	5610125000	3UAX/Lu	200-38422-67	9423 390 20711	P830
Magnesium	Mg	N2025330	5610125100	3QNY/Mg	200-38422-12	9423 390 20121	P831
Manganese	Mn	N2025331	5610125200	3UNX/Mn	200-38422-13	9423 390 20251	P832
Mercury	Hg	N2025322	5610125300	3UNX/Hg	200-38422-28	9423 390 20801	P833
Molybdenum	Mo	N2025332	5610125400	3UNX/Mo	200-38422-31	9423 390 20421	P834
Neodymium	Nd	N2025335	5610125500	3BNX/Nd	200-38422-68	9423 390 20601	P835
Nickel	Ni	N2025336	5610125600	3QNY/Ni	200-38422-15	9423 390 20281	P836
Niobium	Nb	N2025334	5610125700	3UAX/Nb	200-38422-32	9423 390 20411	P837
Osmium	Os	N2025337	5610125800	3UNX/Os	200-38422-69	9423 390 20761	P838
Palladium	Pd	N2025340	5610125900	3UNX/Pd	200-38422-41	9423 390 20461	P839
Platinum	Pt	N2025342	5610126100	3UNX/Pt	200-38422-20	9423 390 20781	P840
Potassium	K	N2025326	5610126200	3BNX/K	200-38422-22	9423 390 20191	P841
Phosphorus	P	N2025338	5610126000	3QNY/P	-	9423 390 20151	P874
Praseodymium	Pr	N2025341	5610126300	3BNX/Pr	200-38422-70	9423 390 20591	P842
Rhenium	Re	N2025344	5610126400	3UAX/Re	200-38422-44	9423 390 20751	P843
Ruthenium	Ru	N2025346	5610126700	3UNX/Ru	200-38422-45	9423 390 20441	P846
Rhodium	Rh	N2025345	5610126500	3UAX/Rh	200-38422-49	9423 390 20451	P844
Rubidium	Rb	N2025343	5610126600	3BNX/Rb	200-38422-33	9423 390 20371	P845
Samarium	Sm	N2025351	5610126800	3UNX/Sm	200-38422-72	9423 390 20621	P847
Scandium	Sc	N2025348	5610126900	3UNX/Sc	200-38422-71	9423 390 20211	P848
Selenium	Se	N2025349	5610127000	3QNY/Se	200-38422-46	9423 390 20341	P849
Silicon	Si	N2025350	5610127100	3QNY/Si	200-38422-16	9423 390 20141	P850
Silver	Ag	N2025300	5610127200	3UAX/Ag	200-38422-02	9423 390 20471	P851
Sodium	Na	N2025333	5610127300	3UNX/Na	200-38422-14	9423 390 20111	P852
Strontium	Sr	N2025353	5610127400	3BAX/Sr	200-38422-34	9423 390 20381	P853
Tantalum	Ta	N2025354	5610127500	3UNX/Ta	200-38422-35	9423 390 20731	P854
Terbium	Tb	N2025355	5610127700	3BNX/Tb	200-38422-73	9423 390 20651	P856
Tellurium	Te	N2025356	5610127600	3QNY/Te	200-38422-47	9423 390 20521	P855
Thallium	Tl	N2025359	5610127800	3UNX/Tl	200-38422-74	9423 390 20811	P857
Thorium	Th	N2025357	-	3UNX/Th	-	9423 390 20901	P858
Thulium	Tm	N2025360	5610128000	3UNX/Tm	200-38422-75	9423 390 20691	P859
Tin	Sn	N2025352	5610128100	3QNY/Sn	200-38422-18	9423 390 20501	P860
Titanium	Ti	N2025358	5610128200	3UNX/Ti	200-38422-17	9423 390 20221	P861
Tungsten	W	N2025362	5610128300	3UNX/W	200-38422-36	9423 390 20741	P862
Vanadium	V	N2025361	5610128500	3UNX/V	200-38422-19	9423 390 20231	P864
Ytterbium	Yb	N2025364	5610128600	3UNX/Yb	200-38422-77	9423 390 20701	P865
Yttrium	Y	N2025363	5610128700	3BNX/Y	200-38422-76	9423 390 20391	P866
Zinc	Zn	N2025365	5610128800	3QNY/Zn	200-38422-23	9423 390 20301	P867
Zirconium	Zr	N2025366	5610128900	3UNX/Zr	200-38422-37	9423 390 20401	P868

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 ©2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.