



7890A/6890 Gas Chromatograph Supplies Quick Reference Guide

Agilent Supplies for Agilent Instruments

Agilent Technologies is committed to optimizing your laboratory's productivity. As part of this long-term commitment, we have designed our supplies to match our instruments precisely. You get consistent results and optimum instrument performance, all day, every day.

Simply mark the supplies that you need or want to keep on hand. Then visit our Web site at www.agilent.com/chem/store for easy and convenient ordering. You can also call (or fax this marked list to) your local Agilent Representative or Agilent Authorized Distributor.

For your convenience, the shopping list below contains the most commonly ordered supplies and parts.

Description	Part No.	Recommended Quantity	Mark for Order
Column Installation Supplies			
Universal column nut	5181-8830	3	<input type="checkbox"/>
Ferrules, 10/pk	Ferrule ID (mm)	Column ID (mm)	
Graphite	0.5	0.1, 0.2, 0.25, 0.32	5080-8853 2 <input type="checkbox"/>
Graphite	1.0	0.53	5080-8773 2 <input type="checkbox"/>
Vespel/Graphite	0.4	0.1, 0.2, 0.25	5181-3323 2 <input type="checkbox"/>
Vespel/Graphite	0.5	0.32	5062-3514 2 <input type="checkbox"/>
Vespel/Graphite	0.8	0.45, 0.53	5062-3512 2 <input type="checkbox"/>
Column ferrule installation tool	19251-80680	1	<input type="checkbox"/>
Diamond-tipped column cutting pencil	420-1000	1	<input type="checkbox"/>
Inlet Septa			
11 mm Non-Stick BTO septa, 50/pk	5183-4757	1	<input type="checkbox"/>
11 mm Advanced Green septa, 50/pk	5183-4759	1	<input type="checkbox"/>
11mm Long Life Inlet septa, 50/pk	5183-4761	1	<input type="checkbox"/>
5 mm through hole septa, 25/pk	5181-1260	1	<input type="checkbox"/>
5 mm Long Life Inlet septa, 50/pk	5183-4762	1	<input type="checkbox"/>
High Pressure Merlin Microseal	5182-3442	1	<input type="checkbox"/>
Merlin Microseal, seal only	5182-3444	1	<input type="checkbox"/>
Split/Splitless Inlet Supplies			
Liner, split, straight, glass wool, non-deactivated	19251-60540	5	<input type="checkbox"/>
Splitless, single taper, deactivated, 4 mm ID	5181-3316	5	<input type="checkbox"/>
Liner, splitless, single-taper, glass wool, deactivated	5062-3587	5	<input type="checkbox"/>
Split, single taper, low pressure drop, glass wool	5183-4647	5	<input type="checkbox"/>
Non-Stick Fluorocarbon O-ring, 10/pk	5188-5365	2	<input type="checkbox"/>
Non-Stick Fluorocarbon Liner O-ring for Flip Top, 10/pk, 6890 only	5188-5366	2	<input type="checkbox"/>
Gold plated seal kit	5188-5367	2	<input type="checkbox"/>

Continued on the next page.



Agilent Analytical Direct 1-800-227-9770 (in the U.S. and Canada)

Description	Part No.	Recommended Quantity	Mark for Order
Cool On-Column Inlet Supplies			
Septum nut base for 320 mm assembly	19245-80521	1	<input type="checkbox"/>
Septum nut base for 530 mm assembly	G1545-80520	1	<input type="checkbox"/>
Spring	19245-60760	1	<input type="checkbox"/>
Inserts			
For 200 µm columns, 1 ring	19245-20510	5	<input type="checkbox"/>
For 250 µm columns, 6 rings	19245-20515	5	<input type="checkbox"/>
For 320 µm columns, 5 rings	19245-20525	5	<input type="checkbox"/>
Inserts for capillary columns	19245-20580	5	<input type="checkbox"/>
Packed Column Inlet Supplies			
O-ring, Fluorocarbon , 12/pk	5080-8898	1	<input type="checkbox"/>
Ferrule, 1/4 in. Vespel	5080-8774	2	<input type="checkbox"/>
1/4 in. nut, brass, 10/pk	5180-4105	1	<input type="checkbox"/>
Disposable glass liner, 170 µl internal volume	5080-8732	1	<input type="checkbox"/>
Flame Ionization Detector (FID) Supplies			
Ignitor glow plug assembly	19231-60680	1	<input type="checkbox"/>
Blanking plug, finger-tight style	5020-8294	1	<input type="checkbox"/>
FID collector assembly	G1531-60690	1	<input type="checkbox"/>
Flapper shaft assembly	G1530-80560	1	<input type="checkbox"/>
High-temperature jet, SimDis 0.018 in. ID tip	G1531-80620	1	<input type="checkbox"/>
Jet, 0.011 in. ID tip, capillary adaptable	19244-80560	1	<input type="checkbox"/>
Jet, packed standard 0.018 in. ID tip	18710-20119	1	<input type="checkbox"/>
FID flow measuring insert	19301-60660	1	<input type="checkbox"/>
FID cleaning kit	9301-0985	1	<input type="checkbox"/>
Nitrogen Phosphorus Detector (NPD) Supplies			
NPD white ceramic bead assembly	G1534-60570	2	<input type="checkbox"/>
NPD black ceramic bead assembly	5183-2007	2	<input type="checkbox"/>
Jet, 0.011 in./0.29 mm ID tip, capillary dedicated	G1531-80560	1	<input type="checkbox"/>
Jet, 0.011 in. ID tip, capillary adaptable	19244-80560	1	<input type="checkbox"/>
FID and NPD Adapters			
FID/NPD adapter for capillary column	19244-80610	1	<input type="checkbox"/>
Ferrule, 1/4 in. Vespel	5080-8774	2	<input type="checkbox"/>
530 µm, 1.0 mm ID graphite ferrule	5080-8773	2	<input type="checkbox"/>
320 µm, 0.5 mm ID graphite ferrule	5080-8853	2	<input type="checkbox"/>
1/4 in. nut, brass	5180-4105	1	<input type="checkbox"/>
Thermal Conductivity Detector (TCD) Supplies			
TCD Capillary Column Adapter	G1532-80540	1	<input type="checkbox"/>
Electron Capture Detector (ECD) Supplies			
Micro ECD wipe test kit	18713-60050	1	<input type="checkbox"/>

Continued on the next page.

Agilent Analytical Direct 1-800-227-9770 (in the U.S. and Canada)

Description	Part No.	Recommended Quantity	Mark for Order
Tools			
GC Installation Kit with Gas Purifiers (no tools)	19199N	1	<input type="checkbox"/>
GC Installation Kit with Tools (no gas purifiers)	19199M	1	<input type="checkbox"/>
General laboratory tool kit	5180-4162	1	<input type="checkbox"/>
Injection Port Cleaning Kit	480-0003	1	<input type="checkbox"/>
Gas Management Supplies			
CGA 580, 125 psig max (8.6 bar), He, Ar, N ₂	5183-4644	1	<input type="checkbox"/>
OT3 Trap	OT3-2	1	<input type="checkbox"/>
Hydrocarbon S-Trap, 1/8 in. fittings	5060-9096	1	<input type="checkbox"/>
Moisture S-trap, preconditioned, 1/8 in. fittings	5060-9084	1	<input type="checkbox"/>
Capillary Flow Technology Supplies			
Ultimate Union kit, deactivated	G3182-61580	1	<input type="checkbox"/>
Internal nut	G2855-20530	1	<input type="checkbox"/>
Swaging nut	G2855-20555	1	<input type="checkbox"/>
Column storage fitting	G2855-20590	2	<input type="checkbox"/>
Fused silica, deactivated, 0.15 mm x 1 m	160-2625-1	1	<input type="checkbox"/>
Fused silica, deactivated, 0.15 mm x 5 m	160-2625-5	1	<input type="checkbox"/>
Fused silica, deactivated, 0.15 mm x 10 m	160-2625-10	1	<input type="checkbox"/>
SilTite Metal Ferrules, 0.20-0.25 mm ID capillary columns	5188-5361	2	<input type="checkbox"/>
SilTite Metal Ferrules, 0.32 mm ID capillary columns	5188-5362	2	<input type="checkbox"/>
SilTite Metal Ferrules, 0.53 mm ID capillary columns	5188-5363	2	<input type="checkbox"/>

Agilent Columns and Supplies – to Help You Get the Results You Need

For greater confidence in your chromatography results, you can optimize your total system with quality Agilent columns and supplies. A wide range of columns, supplies, kits, and accessories is designed, manufactured, and tested to rigorous Agilent specifications, under a quality system registered to ISO 9001. Why risk compromising your analytical results with anything less than genuine Agilent supplies?

For Agilent's complete listing of columns and supplies, visit us online at www.agilent.com/chem/supplies.

Agilent Technologies shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material. Information, description, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2007
Printed in U.S.A. June 4, 2007
5989-6724EN



Agilent Technologies