



7694 Headspace Sampler 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
Vials, Caps and Septa			
Headspace crimp top, flat bottom vials, 20 mL, 100/pk	5182-0837	1	<input type="checkbox"/>
Headspace crimp top, flat bottom vials, 10 mL, 100/pk	5182-0838	1	<input type="checkbox"/>
Silver aluminum crimp caps with 20 mm, PTFE/silicone septa, 100/pk	5183-4477	1	<input type="checkbox"/>
Silver aluminum crimp caps with safety feature, 20 mm, PTFE/silicone septa, 100/pk	5183-4478	1	<input type="checkbox"/>
Silver aluminum crimp caps with 20 mm molded PTFE/butyl septa, 100/pk	5183-4479	1	<input type="checkbox"/>
Silver aluminum crimp caps with safety feature, 20 mm molded PTFE/butyl septa, 100/pk	5183-4480	1	<input type="checkbox"/>
Vial kit: 20 mL Headspace crimp top, flat bottom vials, silver aluminum one-piece crimp caps with safety feature, molded gray PTFE/black butyl septa, 100/pk	5182-0839	1	<input type="checkbox"/>
Vial kit: 20 mL Headspace crimp top, flat bottom vials, silver aluminum one-piece crimp caps with safety feature, PTFE/white silicone septa, 100/pk	5182-0840	1	<input type="checkbox"/>
Needles			
Needle, headspace transfer line, deactivated 0.5 mm OD	2322590004	1	<input type="checkbox"/>
Needle for transfer line, 0.25 mm ID, 0.5 mm OD, nickel	301-016-HSP	1	<input type="checkbox"/>
Needle, headspace transfer line, deactivated 0.7 mm OD	2322590005	1	<input type="checkbox"/>
Needle for transfer line, 0.4 mm ID, 0.8 mm OD, nickel	301-015-HSP	1	<input type="checkbox"/>
Needle assembly vial probe, deactivated	301-220-HSP	1	<input type="checkbox"/>
Needle assembly (vial probe)	301-013-HSP	1	<input type="checkbox"/>
Fittings			
Zero dead volume union	325-045-HSP	1	<input type="checkbox"/>
Bulkhead fitting kit	325-024-HSP	1	<input type="checkbox"/>
Transfer line nut	19258-20830	1	<input type="checkbox"/>
Transfer line ferrule	19258-20870	1	<input type="checkbox"/>
Union FF 6MB, 5-piece set	325-062-HSP	1	<input type="checkbox"/>
Union T6 MB, 5-piece set, brass	325-132-HSP	1	<input type="checkbox"/>
Union T5 MA	325-185-HSP	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
Valves			
Valve, solenoid vent kalrez	3600500001	1	<input type="checkbox"/>
Valve, solenoid vial pressurization	3600500002	1	<input type="checkbox"/>
Pressure regulator	330-086-HSP	1	<input type="checkbox"/>
Flow controller	330-055-HSP	1	<input type="checkbox"/>
Tubing and Transfer Lines			
Sample loop, 1 mL, deactivated	2321700003	1	<input type="checkbox"/>
Sample loop, 1 mL, nickel	321-055-HSP	1	<input type="checkbox"/>
Sample loop, 3 mL, deactivated	2321700004	1	<input type="checkbox"/>
Sample loop, 3 mL, nickel	321-056-HSP	1	<input type="checkbox"/>
Tube, needle, 6-port valve, deactivated	301-212-HSP	1	<input type="checkbox"/>
Tube, needle, 6-port valve, nickel	301-169-HSP	1	<input type="checkbox"/>
Tube, vent-valve stainless steel	301-170-HSP	1	<input type="checkbox"/>
Sensor tube, 125 mm PTFE	321-057-HSP	1	<input type="checkbox"/>
Restrictor, stainless steel	321-002-HSP	1	<input type="checkbox"/>
Transfer line, deactivated, 1 m	301-211-HSP	1	<input type="checkbox"/>
Transfer line, 1 m, nickel	301-152-HSP	1	<input type="checkbox"/>
Transfer line, 80 cm, nickel	301-011-HSP	1	<input type="checkbox"/>
Repair, Leak Test, and OQ/PV Supplies			
Strain relief septum nut	301-205-HSP	1	<input type="checkbox"/>
Oven adaptor for 10 mL vials	301-017-HSP	1	<input type="checkbox"/>
Tray adaptors for 10 mL vials, 25/pk	300-301-HSP	1	<input type="checkbox"/>
Headspace leak test kit	G1888-60701	1	<input type="checkbox"/>
OQ/PV Headspace Sample	5182-9733	1	<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
5989-6723EN



Agilent Technologies