



# Agilent GC Autosampler Supplies (Models 7673/G1513A/7683) 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.

For your convenience, the shopping list below contains the most commonly ordered supplies and parts for the Agilent gas chromatograph autosampler. Simply mark the supplies that you need or want to keep on hand. Then visit our Web site at [www.agilent.com/chem/store](http://www.agilent.com/chem/store) for easy and convenient ordering. You can also call (or fax this marked list to) your local Agilent sales office or authorized Agilent distributor.

Description	Part Number	Recommended Quantity	Mark for Order
<b>Screw-Top Vials and Caps</b>			
Wide opening, clear, screw-top vials (100/pk)	5182-0714	1	<input type="checkbox"/>
Wide opening, amber, screw-top vials, write-on spot (100/pk)	5182-0716	1	<input type="checkbox"/>
Pre-assembled clear vials, blue screw caps, PTFE/silicone septa (100/pk)	5182-0557	1	<input type="checkbox"/>
Non-assembled, 500 vial convenience pack, clear vials, blue screw caps, PTFE/red rubber septa	5182-0867	1	<input type="checkbox"/>
Wide opening, clear, screw-top vials (1000/pk)	5183-2067	1	<input type="checkbox"/>
Blue screw caps, PTFE/red rubber septa	5182-0717	1	<input type="checkbox"/>
Green screw caps, PTFE/red rubber septa	5182-0718	1	<input type="checkbox"/>
Red screw caps, PTFE/red rubber septa	5182-0719	1	<input type="checkbox"/>
Blue screw caps, PTFE/silicone/PTFE septa	5182-0723	1	<input type="checkbox"/>
Green screw caps, PTFE/silicone/PTFE septa	5182-0724	1	<input type="checkbox"/>
Red screw caps, PTFE/silicone/PTFE septa	5182-0725	1	<input type="checkbox"/>
<b>Crimp-Top Vials and Caps</b>			
Wide opening, clear, crimp-top vials (100/pk)	5181-3375	1	<input type="checkbox"/>
Wide opening, amber, crimp-top, write-on spot, 2 ml (100/pk)	5181-3376	1	<input type="checkbox"/>
500 vial convenience pack, clear crimp vials, silver caps, PTFE/red rubber septa	5181-3400	1	<input type="checkbox"/>
Wide opening, clear, crimp-top, 2 ml (1000/pk)	5183-4491	1	<input type="checkbox"/>
Wide opening, clear, crimp-top, write-on spot, 2 ml (1000/pk)	5183-4492	1	<input type="checkbox"/>
Silver aluminum seal, PTFE/red rubber septa (100/pk)	5181-1210	1	<input type="checkbox"/>
Silver aluminum seal, PTFE/silicone septa (100/pk)	5182-0552	1	<input type="checkbox"/>
Silver aluminum seal PTFE/silicone/PTFE septa (100/pk)	5181-1211	1	<input type="checkbox"/>
Blue aluminum seal, PTFE/red rubber septa (100/pk)	5181-1215	1	<input type="checkbox"/>
Green aluminum seal, PTFE/red rubber septa (100/pk)	5181-1216	1	<input type="checkbox"/>
Red aluminum seal, PTFE/red rubber septa (100/pk)	5181-1217	1	<input type="checkbox"/>

Continued on the next page.



Description	Part Number	Recommended Quantity	Mark for Order
<b>Snap-Top Vials and Caps</b>			
Wide opening, clear, snap-top vials (100/pk)	5182-0544	1	<input type="checkbox"/>
Wide opening, amber, snap-top, write-on spot, 2 ml (100/pk)	5182-0545	1	<input type="checkbox"/>
500-vial convenience pack, clear snap vials, polyethylene caps, PTFE/red rubber septa	5182-0547	1	<input type="checkbox"/>
Clear polypropylene snap caps, PTFE septa (100/pk)	5182-0540	1	<input type="checkbox"/>
Clear polypropylene snap caps, PTFE/red rubber septa (100/pk)	5182-0550	1	<input type="checkbox"/>
<b>Polypropylene Vials</b>			
Polypropylene, crimp/snap top, 100 ul (1000/pk)	9301-0978	1	<input type="checkbox"/>
Polypropylene/glass lined, crimp/snap top, 100 ul, (100/pk)	9301-0977	1	<input type="checkbox"/>
Polypropylene, wide opening, crimp/snap top, 1 ml (100/pk)	5182-0567	1	<input type="checkbox"/>
<b>Vial Inserts</b>			
100 uL, mandrel point inserts with polymer feet (100/pk)	5181-1270	1	<input type="checkbox"/>
250 uL, flat-bottom glass inserts (500/pk)	5181-3377	1	<input type="checkbox"/>
<b>Crimping and Decapping Supplies</b>			
Electronic crimper for 2 mL vials with battery and charger	5183-4763	1	<input type="checkbox"/>
Electronic decapper for 2 mL vials with battery and charger	5184-3567	1	<input type="checkbox"/>
Manual Crimp holder without crimp head	5062-2495	1	<input type="checkbox"/>
Crimp head for 12 mm caps (for manual crimp holder)	5001-3755	1	<input type="checkbox"/>
Decapper head for 12 mm caps (for manual crimp holder)	5001-3758	1	<input type="checkbox"/>
<b>Syringes</b>			
Syringe, 10 ul tapered, FN 23-26s/42/HP (6/pk)	5181-3360	1	<input type="checkbox"/>
Syringe, 10 ul straight, FN 23s/42/HP (6/pk)	9301-0725	1	<input type="checkbox"/>
Syringe, 10 ul tapered, FN, gastight, 23-26s/42/HP (6/pk)	5181-3361	1	<input type="checkbox"/>
Syringe, 5 ul tapered, FN 23-26s/42/HP (6/pk)	5181-8810	1	<input type="checkbox"/>
Syringe, 5 ul straight, FN 23s/42/HP (6/pk)	5182-0875	1	<input type="checkbox"/>
Syringe, 5 ul straight, FN 23/42/HP	9301-0892	1	<input type="checkbox"/>
<b>Capillary Inlet Liners</b>			
4 mm id, 900 µl, single taper, deactivated borosilicate, glass wool packing, splitless	5062-3587	2	<input type="checkbox"/>
4 mm id, 900 µl, single taper, deactivated borosilicate splitless	5181-3316	2	<input type="checkbox"/>
4 mm id, 990 µl, split, borosilicate, glass wool packing	19251-60540	2	<input type="checkbox"/>
Split, low pressure drop, taper, glass wool, deactivated	5183-4647	2	<input type="checkbox"/>
General purpose, split/splitless, taper, glass wool, deactivated	5183-4711	2	<input type="checkbox"/>

Continued on the next page.

**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 for the Agilent GC autosampler. 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 consumables?

For Agilent's complete listing of columns and supplies, visit us on the World Wide Web at [www.agilent.com/chem/supplies](http://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. 2005

Printed in U.S.A. April 6, 2005  
5989-2196EN



**Agilent Technologies**