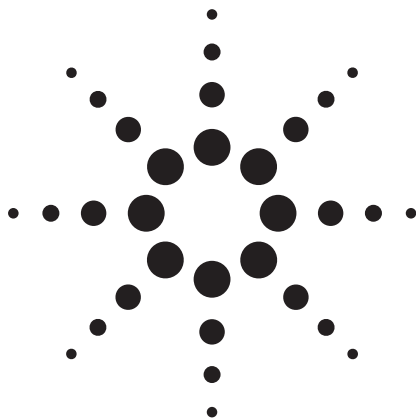




Agilent 5975/5973 MSD Supplies

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.

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

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 sales office or authorized Agilent distributor.

Description	Part No.	Recommended Quantity	Mark for Order
Cleaning and Maintenance Supplies			
One Year Maintenance Kit (for diffusion pump systems)	5183-2096	1	☐
MSD Tool Kit, 5975/5973	G1099-60566	1	☐
Nylon gloves, lint-free, large	8650-0030	4	☐
Abrasive sheets, 5/pk	5061-5896	2	☐
Cloths, lint free, 15/pk	05980-60051	2	☐
Cotton swabs, 100/pk	5080-5400	2	☐
Santovac 5P Ultra diffusion pump fluid, 18.5 ml	6040-0809	2	☐
Foreline pump oil, P3, 0.5 liter	6040-0621	1	☐
Ion Source Supplies			
Filament, EI High Temperature for 5975	G2590-60053	2	☐
Filament, 5975/5973 CI	G1099-80053	1	☐
Lens insulator, 2/pk	05971-20130	1	☐
MS interface column nut	05988-20066	1	☐
Repeller Assembly Supplies			
Repeller insulator	G1099-20133	2	☐
Inlet Septa			
11 mm Non-Stick BTO septa, 50/pk	5183-4757	2	☐
Merlin Microseal			
High pressure Merlin Microseal starter kitIncludes microseal septum and nut	5182-3442	2	☐
Microseal high pressure septum	5182-3444	1	☐
Ferrules			
0.4 mm Vespel Graphite ferrule, for 200/250 µm columns, 10/pk	5062-3508	1	☐
0.5 mm Vespel Graphite ferrule, for 320 µm columns, 10/pk	5062-3506	1	☐
250 µm Vespel/Graphite ferrule, 10/pk	5181-3323	1	☐
SilTite Metal Ferrules, 1/16 in.Includes 2 column nuts	5184-3571	1	☐
SilTite Metal Ferrules, 1/16 in. x 0.4mm IDIncludes 2 column nuts	5184-3569	1	☐
SilTite Metal Ferrules, 1/16 in. x 0.5mm IDIncludes 2 column nuts	5184-3570	1	☐

Continued on the next page.



Agilent Technologies

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

Description	Part No.	Recommended Quantity	Mark for Order
Detector Supplies			
Electron multiplier replacement horn	05971-80103	1	☐
Miscellaneous Supplies			
Gas leak detector, 115V	5182-9646	1	☐
ADM 1000	220-1170	1	☐
Big Universal Trap, 1/8 in. fittings	RMSH-2	1	☐
PFTBA MS Sample Kit, 0.5 ml	05971-60571	1	☐
Replacement glass bulb for PFTBA test sample (5975)	G3170-80002	1	☐
PFDTD sample	8500-8130	1	☐
Replacement glass vial for PFTBA test sample (5973)	05980-20018	1	☐

Description	Non-Inert Part No.	Mark for Order	Inert Part No.	Mark for Order
5975/5973 MSD Ion Source Parts (EI)				
Ion source body	G1099-20130	☐	G2589-20043	☐
Repeller	G1099-20132	☐	G2589-20044	☐
Interface socket	G1099-20136	☐	G1099-20136	☐
Drawout plate, 3 mm	05971-20134	☐	G2589-20100	☐
Drawout plate, 6 mm	G3163-20530	☐	G2589-20045	☐
Drawout cylinder	G1072-20008	☐	G1072-20008	☐
Ion focus lens	05971-20143	☐	05971-20143	☐
Entrance lens	05971-20126	☐	G3170-20126	☐
Repeller insulator	G1099-20133	☐	G1099-20133	☐
Set screw for lens stack	0515-1446	☐	G1999-20022	☐
Washer, for Repeller M3	3050-0891	☐	3050-0891	☐
Nut, 5.5 mm	0535-0071	☐	0535-0071	☐
Ion source sensor	G1099-60104	☐	G1099-60104	☐
Lens insulator, 2/pk	05971-20130	☐	05971-20130	☐
Ion source assembly	G1099-60106	☐	G3170-65710	☐
Anodized source heater assembly	G3169-60177	☐	G3169-60177	☐
Repeller assembly	G1099-60170	☐	G2589-60102	☐
Screw for filament on the source	0515-1046	☐	G1999-20021	☐

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 on the World Wide Web 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
5989-5383EN



Agilent Technologies