



# Agilent Inert Sample Introduction Kit for HF Digests with Axial ICP-OES

Part number 9910126000

## Contents List

Description	Quantity	Part Number
Barb connector for waste line, 1/16 inch to 1/16 inch	5	1610093000
Barb connector for waste line, 1/16 inch to 1/8 inch	5	1610088600
Capillary tubing for sample inlet, connects to peristaltic pump tubing	2 m	2410020500
Tubing, spray chamber to peri-pump waste line, suits Sturman-Masters spray chamber	1 m	3710033900
Tubing, spray chamber to peri-pump waste line, suits cyclonic spray chamber	1 m	3710024600
Transfer tubing, spray chamber to torch	1 m	3710033400
Connecting tubing, to join pump tubes to capillary tubing	0.5 m	3710033700
Peristaltic pump tubes, PVC, blue/blue, 12/pk	1	3710034600
Peristaltic pump tubes, PVC, white/white, 12/pk	1	3710034400
Bracket for mounting Sturman-Masters spray chamber on axial ICP	1	410328600
V-groove nebulizer assembly	1	9910057400
Instruction sheet, V-groove nebulizer	1	
Nebulizer cleaning wire for V-groove nebulizer, 0.15 mm diameter	0.3 m	
Sturman-Masters spray chamber assembly, Ertalyte	1	110593190
Holder nebulizer, Ertalyte	1	
Semi-demountable torch with 1.8 mm id alumina injector	1	9910084700

Agilent leads the way in spectroscopy innovation. Our comprehensive and trusted portfolio offers you the most diverse application coverage for AA, ICP-OES, and ICP-MS, while our unique MP-AES and triple quadrupole ICP-MS technologies deliver new possibilities for your laboratory.

Maximize your productivity and data quality with genuine Agilent parts and supplies. Visit the Agilent website at [www.agilent.com/chem/ICP-OESsupplies](http://www.agilent.com/chem/ICP-OESsupplies) to find the recommended supplies for your instrument.



**Agilent Technologies**

[www.agilent.com/chem/ICP-OESsupplies](http://www.agilent.com/chem/ICP-OESsupplies)

Agilent 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, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2014  
Printed in Malaysia  
August 7, 2014  
5991-5098EN



**Agilent Technologies**