



# Agilent Sample Introduction Kit for Common Organic Solvents with Axial ICP-OES

Part number 9910126300

## 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
One-piece low-flow torch for organic solvents with axial ICP, 1.4 mm id injector	1	2010104800
Conikal glass concentric nebulizer	1	2010106800
Capillary tubing for sample inlet, connects to peristaltic pump tubing	2 m	2410020500
Peristaltic pump tubes, PVC Solvaflex, black/black, 12/pk	1	3710034800
Peristaltic pump tubes, PVC Solvaflex, white/white, 12/pk	1	3710035000
Tubing, spray chamber to peri-pump waste line, solvent resistant	1 m	3710035400
Double-pass, glass cyclonic spray chamber, suitable for all ICP-OES	1	7910043700
Bracket for mounting glass cyclonic spray chamber on axial ICP	1	110665200
Transfer tubing, FEP, glass cyclonic spray chamber for axial ICP	1	7910051590
Sample inlet, EzyFit fitting, for concentric glass nebulizer on axial ICP-OES, 0.75 mm id x 1.3 mm od capillary tubing, 10/pk	1	9910107700

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  
5991-5101EN



**Agilent Technologies**