



# Agilent Organic Solvent Tubing Kit for Radial ICP-OES with Inert Sample Introduction System

Suitable for use with low volatile organic solvents

Part number 9910124500

## Contents List

Description	Quantity	Part Number
Barb connector for waste line, 1/16 inch to 1/8 inch	5	1610088600
Barb connector for waste line, 1/16 inch to 1/16 inch	5	1610093000
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
Solvent resistant tubing, spray chamber to peri-pump waste line, suits inert Sturman-Masters spray chamber	1 m	3710035300
Solvent resistant tubing, spray chamber to peri-pump waste line, suits glass cyclonic spray chamber	1 m	3710035400
Transfer tube, FEP, Sturman-Masters spray chamber for radial ICP	1	7910051490

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/specsuppliesinfo](http://www.agilent.com/chem/specsuppliesinfo)

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  
5991-5093EN



**Agilent Technologies**