



Agilent Aqueous Tubing Kit for Axial ICP-OES with Concentric Nebulizer and Glass Cyclonic Spray Chamber

Part number 9910123600

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 glass cyclonic spray chamber	2 m	3710024600
Transfer tubing, FEP, cyclonic spray chamber for axial ICP	1	7910051590
Peristaltic pump tubes, PVC, blue/blue, 12/pk	1	3710034600
Peristaltic pump tubes, PVC, white/white, 12/pk	1	3710034400
Connecting tubing, to join pump tubes to capillary tubing	0.5	3710033700
Sample inlet, EzyFit, for concentric glass nebulizer on axial ICP-OES, 0.75 mm id, 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 to find the recommended supplies for your instrument.



Agilent Technologies

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
5991-5086EN

