



Agilent Aqueous Tubing Kit for Radial ICP-OES with Inert Sample Introduction System

Part number 9910123900

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 glass cyclonic spray chamber	1 m	3710024600
Transfer tubing, FEP, Sturman-Masters spray chamber for radial ICP	1	7910051490
Peristaltic pump tubes, PVC, blue/blue, 12/pk	1	3710034600
Peristaltic pump tubes, PVC, gray/gray, 12/pk	1	3710034500
Connecting tubing, to join pump tubes to capillary tubing	0.5 m	3710033700

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
5991-5089EN



Agilent Technologies