



Agilent Sample Compartment and Torch Box Tubing Kit for ICP-OES

Part number 9910057100

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, to connect to peristaltic pump tubing	2 m	2410020500
Gas supply tubing, to connect argon gas supply from outlet on instrument to plasma/auxiliary gas inlets on torch	1 m	2410023800
Tubing, spray chamber to peri-pump waste line, suits glass cyclonic spray chamber	2 m	3710024600
Tubing, spray chamber to peri-pump waste line, suits Sturman-Masters spray chamber	1 m	3710033900
Tubing, solvent resistant, spray chamber to peri-pump waste line, suits glass cyclonic spray chamber	1.5 m	3710035400
Transfer tubing, spray chamber to torch	0.5 m	3710033400
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
5991-5083EN



Agilent Technologies