



Agilent Volatile Organic Solvents Sample Introduction Kit for Axial ICP-OES

Part number 9810039690

Contents List

Description	Quantity	Part Number
Double-pass glass cyclonic spray chamber	1	7910043700
Bracket for mounting glass cyclonic spray chamber on axial ICP	1	110665200
SeaSpray concentric glass nebulizer, includes EzyFit sample connector, 0.75 mm id × 700 mm, and EzyLok gas connector	1	2010096400
Nebulizer gas connector with connecting tubing and 2 tubing clamps	1	
Peristaltic pump tubes, PVC, white/white, 12/pk	1	3710034400
Transfer tube, FEP, cyclonic spray chamber for axial ICP	1	7910051590

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
5991-5081EN



Agilent Technologies