



# Agilent High Solids Sample Introduction Kit for Radial ICP-OES

Part number 9810039590

## Contents List

Description	Quantity	Part Number
Double-pass glass cyclonic spray chamber	1	7910043700
Bracket for mounting glass cyclonic spray chamber on radial ICP	1	110611700
SeaSpray concentric glass nebulizer, with 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 radial ICP	1	7910051290

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  
5991-5080EN



**Agilent Technologies**