



# Agilent Inert Hi-Solids Sample Introduction Kit for Radial ICP-OES

Part number 9810039490

## Contents List

| Description  | Quantity | Part Number |
|--|----------|-------------|
| Sturman-Masters spray chamber assembly - Ertalyte                  | 1        | 110593190   |
| Holder Nebulizer - Ertalyte  | 1        | 810128100   |
| Bracket for mounting Sturman-Masters spray chamber on radial ICP   | 1        | 410298400   |
| V-groove nebulizer assembly  | 1        | 9910057400  |
| Instruction Sheet V-groove Nebulizer                               | 1        |             |
| Nebulizer cleaning wire for V-groove nebulizer 0.15 mm diameter    | 3        |             |
| Nebulizer gas connector with connecting tubing and 2 tubing clamps | 1        |             |
| Peristaltic pump tubes, PVC, gray/gray. 12/pk                      | 1        | 3710034500  |
| Transfer tube FEP Sturman-Masters spray chamber for radial ICP     | 1        | 7910051490  |

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/ICP-OESsupplies](http://www.agilent.com/chem/ICP-OESsupplies)

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 the USA  
July 22, 2014  
9810039490

