



Agilent Inert Sample Introduction Kit for HF Digests with 5100 Radial View ICP-OES

Part number G8010-68006

Contents List

Description	Quantity	Part Number
Easy-fit, inert demountable torch for 5100 RV ICP-OES, for HF digests, with 1.8 mm id alumina injector and replaceable outer/intermediate tubes. Suitable for radial viewing	1	G8010-60234
Inert double-pass spray chamber with Helix nebulizer seal, ball joint socket, and UniFit drain outlet	1	G8010-80034
Inert OneNeb concentric nebulizer with sample connector and EzyLok gas connector	1	2010126900
Peristaltic pump tubes, PVC Solvaflex, black/black, 12/pk	1	3710034800
Peristaltic pump tubes, PVC Solvaflex, gray/gray, 12/pk	1	3710035200
Clamp for ball joint connector between spray chamber and inert torch, 1/pk	1	G8010-60227

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/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 Malaysia
August 7, 2014
5991-5096EN



Agilent Technologies