



Agilent Sample Introduction Kit for Volatile Organic Solvents with 5100 Dual View ICP-OES

Part number G8010-68005

Contents List

Description	Quantity	Part Number
Easy-fit, demountable torch for 5100 DV ICP-OES, for volatile organics, with 0.8 mm id injector and replaceable outer/intermediate tubes. Suitable for axial and radial viewing	1	G8010-60235
Conikal glass concentric nebulizer, U-series, with UniFit sample connector and Ezylok gas connector	1	G8010-60270
Peristaltic pump tubes, Viton, white/white, 12/pk	1	3710043700
Peristaltic pump tubes, Viton, blue/blue, 12/pk	1	3710043900
Tubing, spray chamber to peri-pump waste line, solvent resistant	2 m	3710035400

Requires, but does NOT include a cooled spray chamber. The IsoMist Peltier-cooled spray chamber is recommended.

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
5991-5095EN



Agilent Technologies