



Agilent Internal Standard Kit Number 1 for ICP-OES

Part number 9910124000

Contents List

Description	Quantity	Part Number
Y-barb connector, 1/16 inch, to tee the internal standard line to the sample line	5	1610132400
Connecting tubing for sample inlet, to join pump tubes to capillary tubing	0.1 m	3710033700
Peristaltic pump tubes, PVC, black/black, 12/pk	1	3710027200
Sample inlet, EzyFit, for concentric glass nebulizer on axial ICP-OES, 0.75 mm id × 1.3 mm od capillary tubing, 10/pk	1	9910107700

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
5991-5090EN

