



# Agilent Sample Compartment O-ring Kit for ICP-OES

Part number 9910057200

## Contents List

Description	Quantity	Part Number
O-ring for nebulizer inlet of glass cyclonic spray chamber, nitrile rubber, 1/4 inch id, 3/8 inch od, 1/16 inch thick	2	6910001200
O-ring for drain and outlet fittings on Sturman-Masters spray chamber, nitrile rubber, 11/16 inch id, 13/16 inch od, 1/16 inch thick	4	6910006000
O-ring for nebulizer holder of the Sturman-Masters spray chamber, nitrile rubber, 7/32 inch id, 11/32 inch od, 1/16 inch thick	2	6910008200
O-ring for top to bottom seal of the Sturman-Masters spray chamber, nitrile rubber, 2-3/16 inch id, 2-3/8 inch od, 3/32 inch thick	2	6910017400
O-ring for nebulizer inlet of the Sturman-Masters spray chamber, nitrile rubber, 7/16 inch id, 5/8 inch od, 3/32 inch thick	2	6910019800
O-ring for injector holder of the semi-demountable torch, nitrile rubber, 5/32 inch id, 9/32 inch od, 1/16 inch thick	2	6910006700

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  
5991-5084EN



**Agilent Technologies**