



Red Pass Filter Replacement Kit For NOA Models 270B and 280

Required Tools

- Small Phillips Head Screwdriver or
- Allen Wrench for Reaction Cell PMT Cap
- Water or alcohol

Procedure

In conjunction with the Reaction Cell cleaning instructions in the Agilent 355 Sulfur/255 Nitrogen Chemiluminescence Detectors - Operation and Maintenance manual, use this kit to replace the IR Pass Filter and O-Rings.

For more information on our products and services, visit our Web site at www.agilent.com/chem.



Copyright © 2007
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610
Printed in the USA May 2007



G6600-90036