



UV Pass Filter (350B) Replacement Kit

Accessory G6600-67004

TOOLS NEEDED

- Small Phillips head screwdriver/Allen Wrench
- Water or similar solvent

QTY	DESCRIPTION	PART #
1	UV PASS FILTER, 1.668" dia x 3MM THICK, 350B	G6600-80032
1	O-RING, 1.3614" ID	G6600-80050
1	O-RING, 1.301" ID	G6600-80051
1	(this procedure)	G6600-90037

In conjunction with the instructions in the Service Manual on Reaction Cell Cleaning, use this kit to replace the UV 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 Feb 2007



G6600-90037