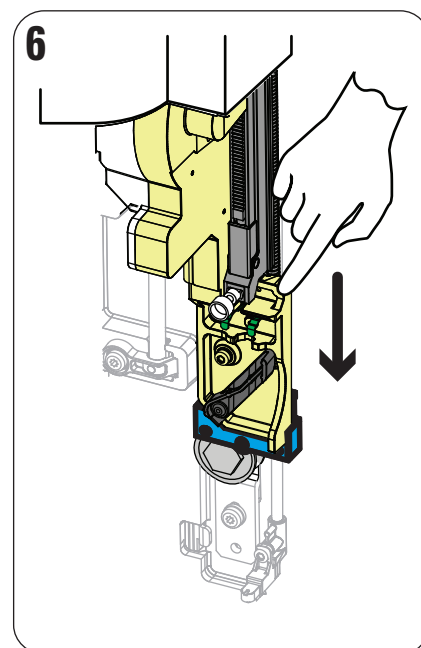
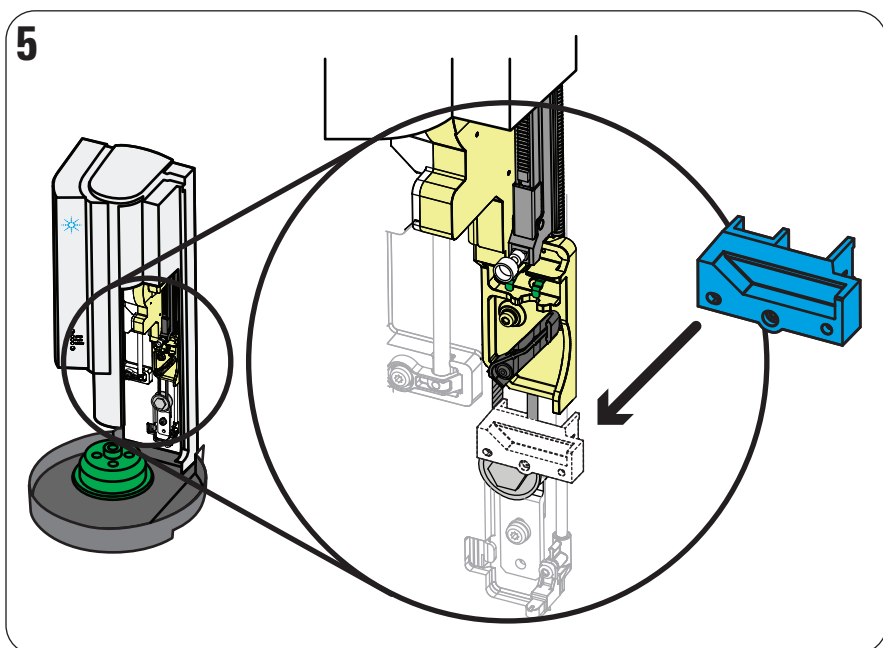
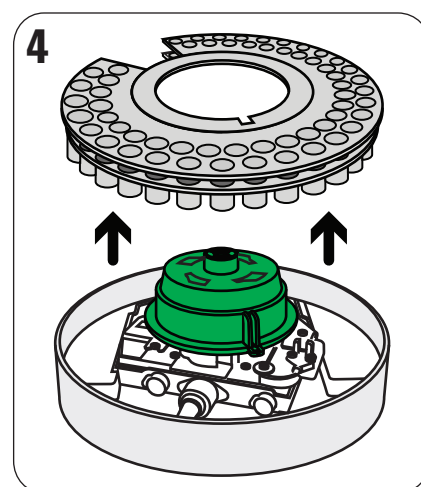
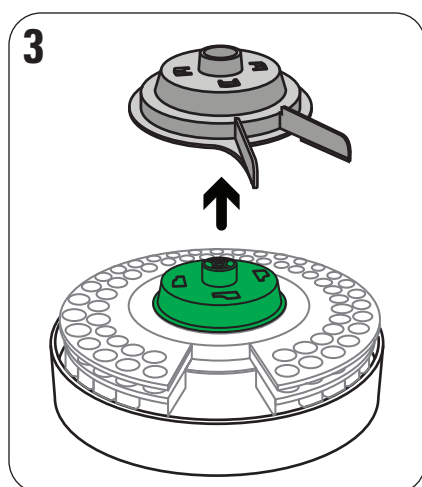
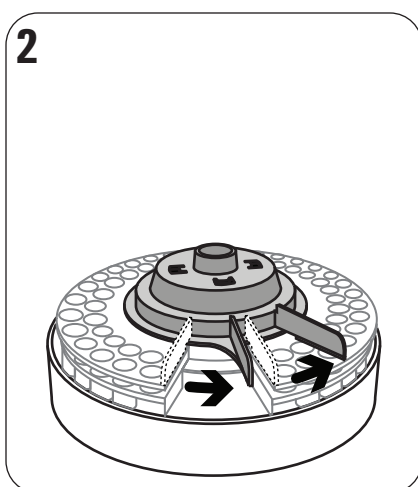
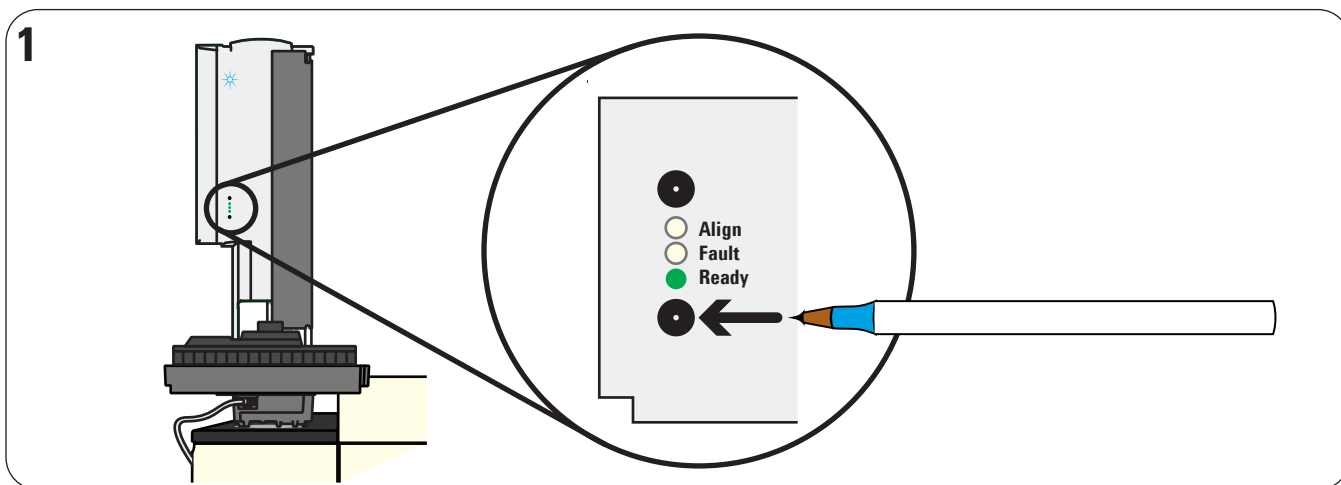
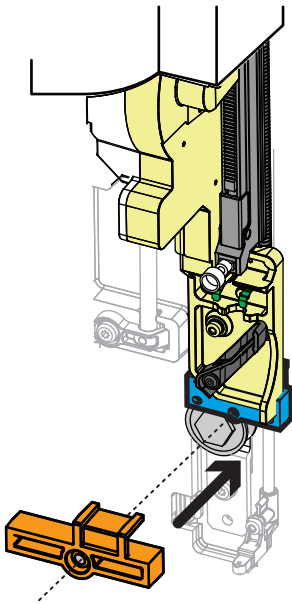
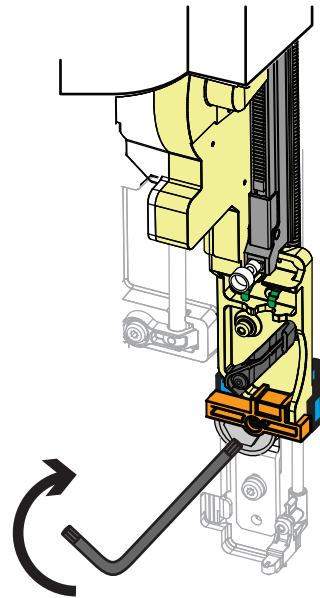




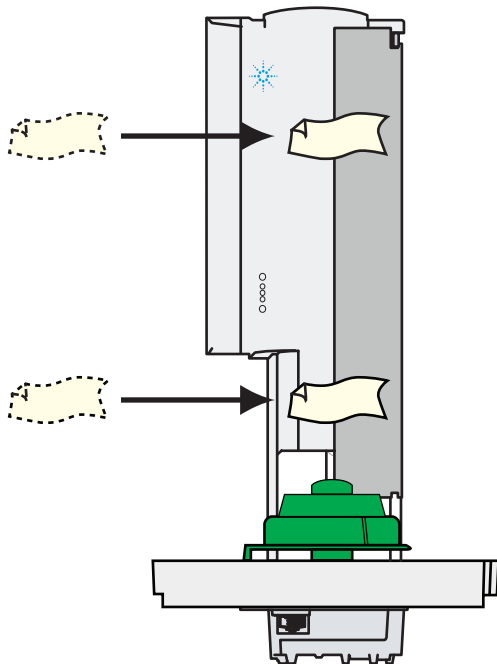
7



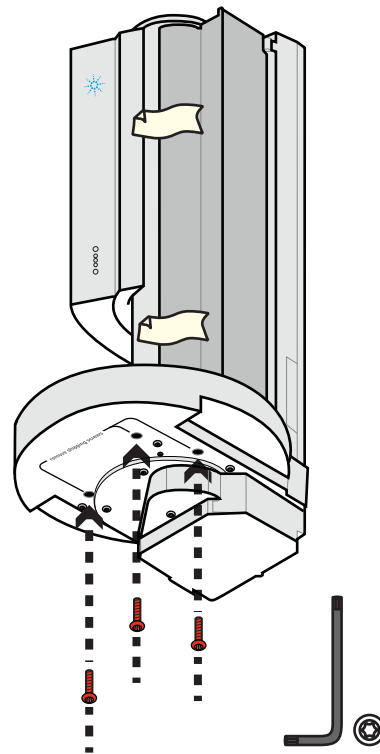
8



9



10



© Agilent Technologies, Inc.
Printed in China, August 2012



G4567-90500 REV A



Agilent Technologies