

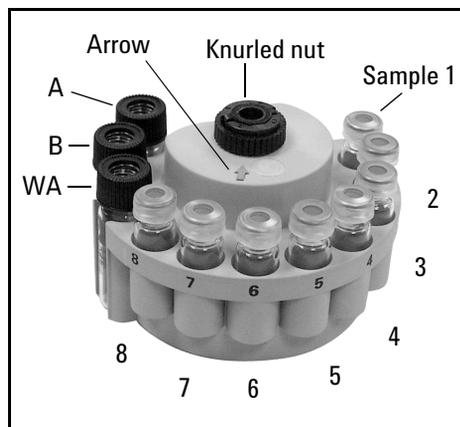
Installing the 8-Vial Turret for the Automatic Liquid Sampler

Agilent 7673A Automatic Liquid Sampler

Accessory G2306A

Introduction

You have just received the G2306A eight-vial turret used on the automatic liquid sampler. This eight-vial sample turret has two solvent positions (labeled A and B), and one waste position (labeled WA). You can control all necessary parameters using either the 6890 series GC keyboard, the Multitechnique ChemStation, the MSD ChemStation, the G1513A injector module rotary switch, or the 3396 integrator. You cannot, however, use the 18596 100-vial sampler tray with this turret.



Hardware Requirements

- G1513A Injector.
- G1512A Controller, firmware A.01.06 or greater.
- Serial number > 3509A01280 or > 3520G00625.
- Tray disabled or not installed.



Agilent Technologies

Software Requirements

- G2070AA, ChemStation, Rev 4.01 or later.
- G1701AA, MSD ChemStation, Rev A.02.00 or later.

Installing the Turret

Use the following instructions to install your new 8-vial turret:

1. Turn off the controller if the injector is installed.
2. Open the tower door. Unscrew and remove the knurled nut from the top of the turret.
3. Rotate the turret so its open section faces the back of the tower, and so the arrow on top of the turret points directly back. Push the stripper arm toward the back, lift the turret up so it clears the center shaft, then remove it from the front of the tower.

Save your standard turret for future use with the 100-vial sampler tray. The 8-vial turret is *incompatible* with the tray.

4. To install the new turret, align the key on the motor shaft with the slot located on the underside of the turret.
5. Push the stripper arm to the back of the tower, insert the turret with its open section facing the back of the tower and arrow pointing back, and install the turret with the key inserted into the slot. Seat the turret onto the shaft, rotating it gently if necessary, until it drops completely into position. The top of the threaded shaft should be almost level with the top of the turret, and the stripper arm should clear the top by approximately 1 mm as the turret is rotated.
6. Replace the knurled nut and tighten firmly (finger-tight).
7. Restore power and check for the green tower ready light. Note that the align LED is on. *Before you start a run* you must perform the syringe carriage height calibration procedure to ensure proper needle depth into the vial. Your instrument *will not run* without this calibration. This procedure (see following) must be performed each time the turret is changed.

Alignment procedure

Each time you change a turret, it is necessary to perform the following alignment procedure:

1. Turn the controller off.
2. Remove the syringe if installed.
3. Remove the needle support rod. If you need help, refer to your *Automatic Liquid Sampler Operating Manual*.
4. Enter diagnostic mode by simultaneously pressing START and STOP buttons while turning on the controller. FAULT, RUN, and READY LEDs (red, yellow, and green respectively) will start lighting in rotation, indicating you are in diagnostic mode.
5. Press the STOP button once. The RED LED should start flashing twice indicating you are in syringe carriage test mode.
6. Press the STOP button three more times until the RED LED is flashing five times per cycle. This indicates you are in alignment mode.
7. Press the START button to begin the alignment procedure. The injector will go through the following steps:
 - The turret will home.
 - The syringe carriage will move up to home, then will move down 240 motor steps. This will be repeated several times with a motor step added each time it moves down until it hits the hard stop on the base assembly. This determines a more precise distance between the syringe carriage and base of the injector. This could take up to 1 minute.
 - The syringe carriage will move up until slightly above the turret, and the turret will move over the inlet.
 - The syringe carriage will move down a step at a time until it hits the top of the turret. This determines a more precise distance between the syringe carriage and top of the turret.
 - The RED LED will return to the beginning of the diagnostic set and flash twice.
8. Replace the needle support rod.

9. Power cycle the controller to exit diagnostic mode and return to the ready state.

Copyright© 2000
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610
Printed in USA JUL 2000



G2306-90107