

Exchanging your Injector

G4567A

| | |
|-------------------------------------|---|
| Safety Instructions | 1 |
| Uninstalling the defective injector | 2 |
| Installing the replacement injector | 2 |
| General | 3 |

Safety Instructions

Always follow your laboratory's safety regulations, and the warnings given in the module's manual.

CAUTION

Condensation within the module will harm the system electronics.

- ✓ Do not store, ship or use your module under conditions where temperature fluctuations could cause condensation within the module.
- ✓ If your system was shipped in cold weather, leave it in its box and allow it to warm slowly to room temperature to avoid condensation.

WARNING

Toxic, flammable and hazardous solvents and reagents

The handling of solvents and reagents can hold health and safety risks.

- ✓ When working with these substances, respect appropriate safety procedures (for example, by wearing goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the vendor, and follow good laboratory practice.

Uninstalling the defective injector

- 1 Turn off the GC.
- 2 Remove the **syringe** and all **samples, wash bottle** and **waste bottle** from the injector turret.
- 3 Unplug the injector power cable from the injector.
- 4 Lift the Injector from the GC and put it aside.

Installing the replacement injector

The injector is shipped with the Std needle-guide foot installed. The CoC foot is shipped with the injector but not installed.

- 1 Unpack the Instrument Exchange Program (IEP) injector from the shipping box.

NOTE

Do not discard any of the packaging; this will be used to return the defective unit to Agilent.

- 2 Remove the three red shipping screws from the bottom of the injector and put them aside.

NOTE

It is important that you remove the shipping screws but keep them for later reuse.

- 3 Remove the shipping tape from injector door.
- 4 Remove the orange/blue shipping clamp from the syringe carriage with a Torx-10 tool.
- 5 Verify that the configuration of the replacement injector is correct, for example, syringe carriage (**Std or LVI**).
- 6 **Cool-on-Column:** (If you do not use CoC, than skip this step)
 - a Use a Torx-10 tool to remove the Std foot.
 - b Remove the T-10 screw and install the CoC foot in the same position.
- 7 Install the injector on the GC post.
- 8 Connect the injector power cable to the injector.
- 9 Turn the GC on and check if the **green light** on the Injector is on.

NOTE

If the injector starts with a red or flushing light, check the following:

- 1 **Red light:** Is the door closed?
- 2 **Red and yellow lights:** Is the injector sitting correctly on the GC post?
- 3 If this does not help, start the alignment procedure by pushing the align button (**ensure that no syringe or bottles are installed before you start**).

General

- 1 Install the **orange/blue shipping clamp** and the **three red shipping screws** from the replacement injector in the defective unit.
- 2 Using the original packing material, pack the defective unit into the shipping box from the exchange unit.
- 3 Fill out the attached "Environmental Health and Safty Sheet (**EHS**)" and place it in the shipping box.
- 4 **Return** the defective unit to Agilent **within five days**.



Discover Agilent