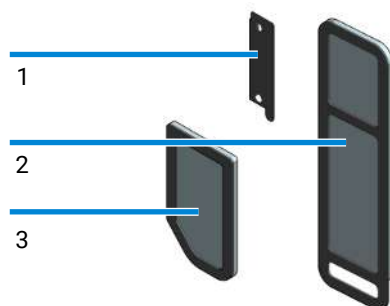


Agilent InfinityLab
 Dust Filter Kit
Installation

This Technical Note describes the installation of the 5720-0026 (Dust filter kit) in an Infinity II/III autosampler.

Overview of Components



5720-0026 (Dust filter kit) consists of:

#	p/n	Description
1		Magnetic cover for electronic box
2		Magnetic filter pad, left
3		Magnetic filter pad, right

CAUTION

Magnetic component

NOTE

The 5720-0026 (Dust filter kit) is compatible with all Agilent Infinity II/III autosamplers.

NOTE

Agilent recommends a yearly replacement of the dust filters.

NOTE

While installing dust filters can help prevent future exposure, it is important to note that they cannot eliminate the risk associated with existing dust contamination already present in the system.



Install the Dust Filter Kit

Install the Dust Filter Kit


When

- The autosampler is exposed to significant amounts of dust or other airborne particles. Dust or other airborne particles can potentially reduce the lifespan of the electronic boards.

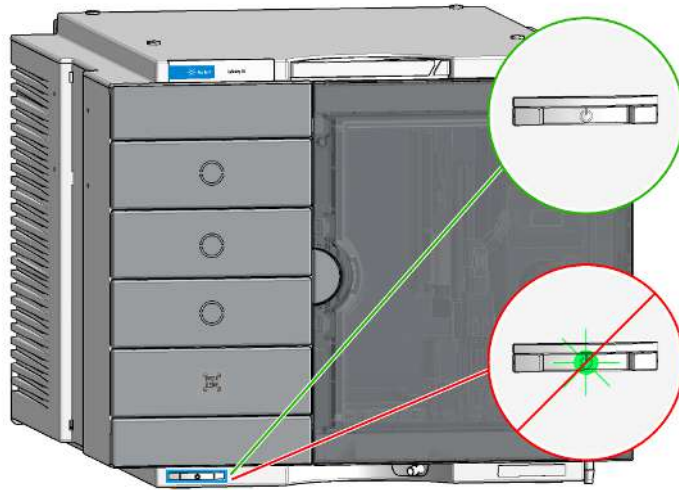
Tools required

Qty.	p/n	Description
1	 5182-3466	Torx screwdriver T10
1	 5023-3089	Torx key set (Part of the G7120-68708 InfinityLab System Tool kit)

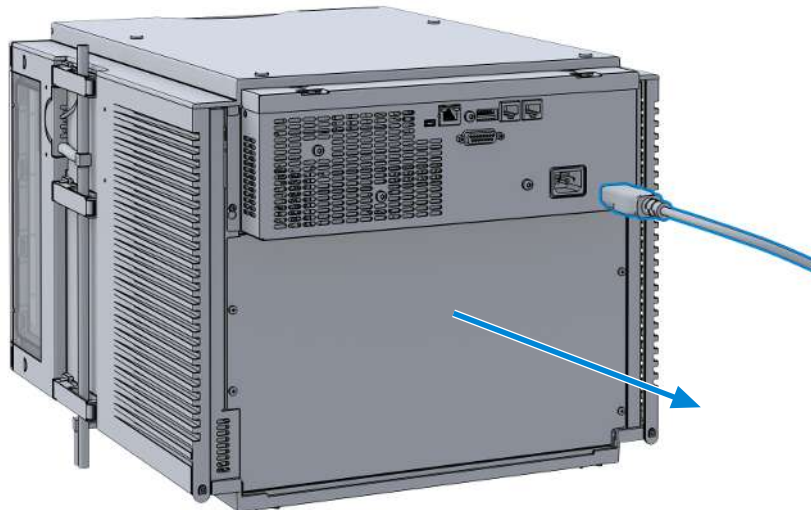
Parts required

Qty.	p/n	Description
1	 5720-0026	Dust filter kit

1 Ensure that the on/off switch on the front of the module is OFF (switch stands out).



- 2 Disconnect the power cable from the sampler.

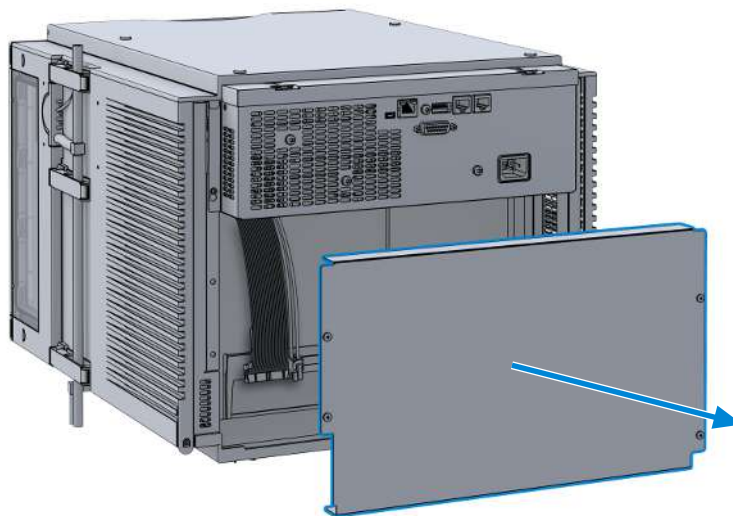


Install the Dust Filter Kit

- 3 Loosen the four screws on the rear of the module.
NOTE: If a Sample Thermostat is installed in the sampler, remove it by following the procedure outlined in the module-specific user manual. Then, continue with step 5.

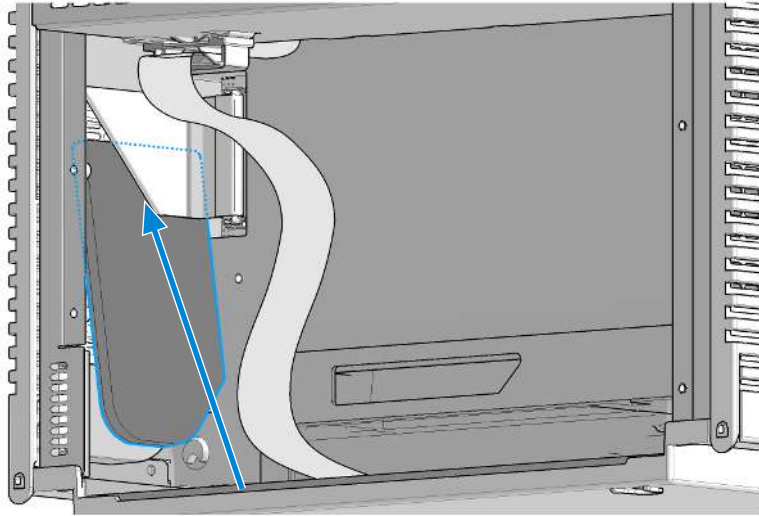


- 4 Remove the sheet metal back cover of the sampler.

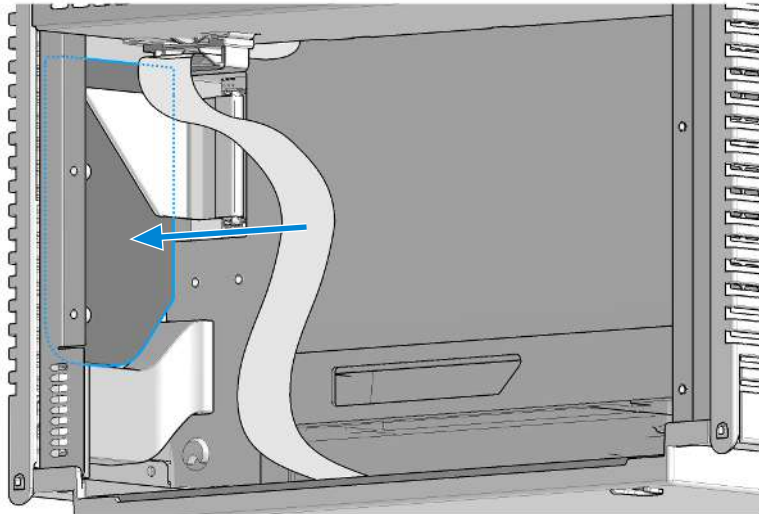


Install the Dust Filter Kit

- 5 Route the upper edge of the smaller magnetic filter pad along the inner surface of the sheet metal housing where the thermostat cables are located until resistance is felt. This area is on the left side when viewed from the back of the module.

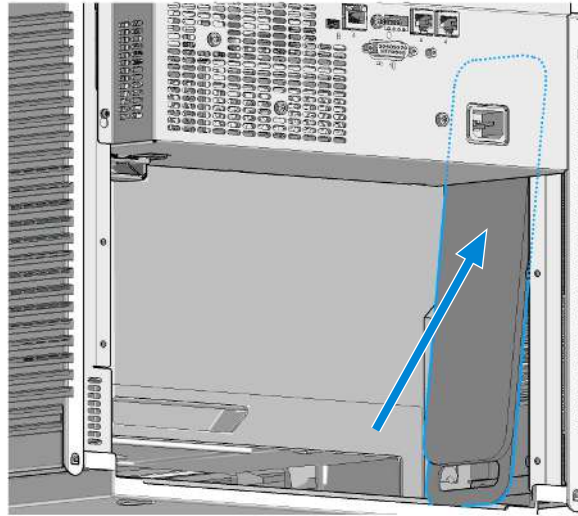


- 6 Adhere the filter pad to the surface, ensuring it is oriented correctly and fully covers all venting holes in the sheet metal housing.

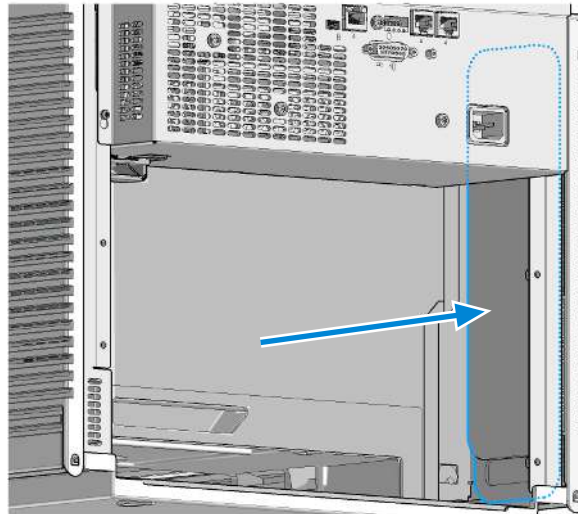


Install the Dust Filter Kit

- 7 Route the upper edge of the larger magnetic filter pad along the inner surface of the sheet metal housing where the flat flexible cable is located until resistance is felt. This area is on the right side when viewed from the back of the module.

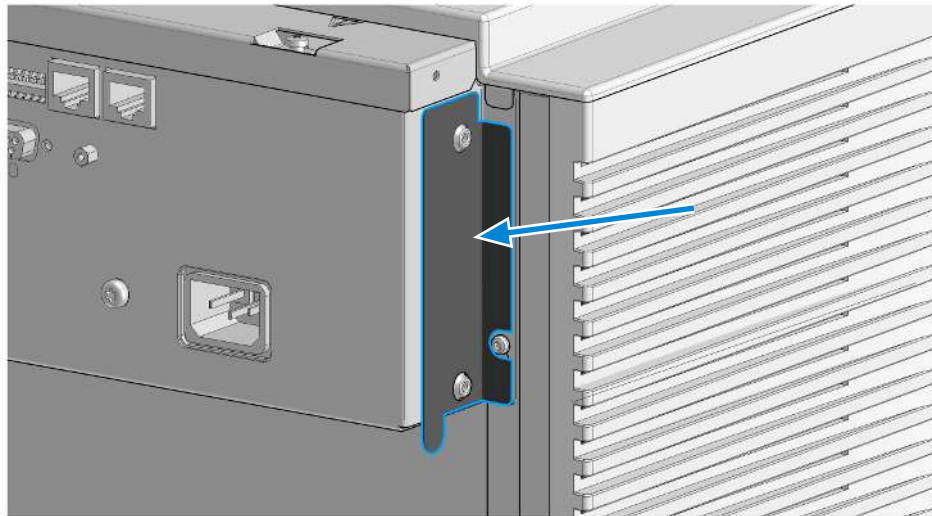


- 8 Adhere the filter pad to the surface, ensuring it is oriented correctly and fully covers all venting holes in the sheet metal housing.

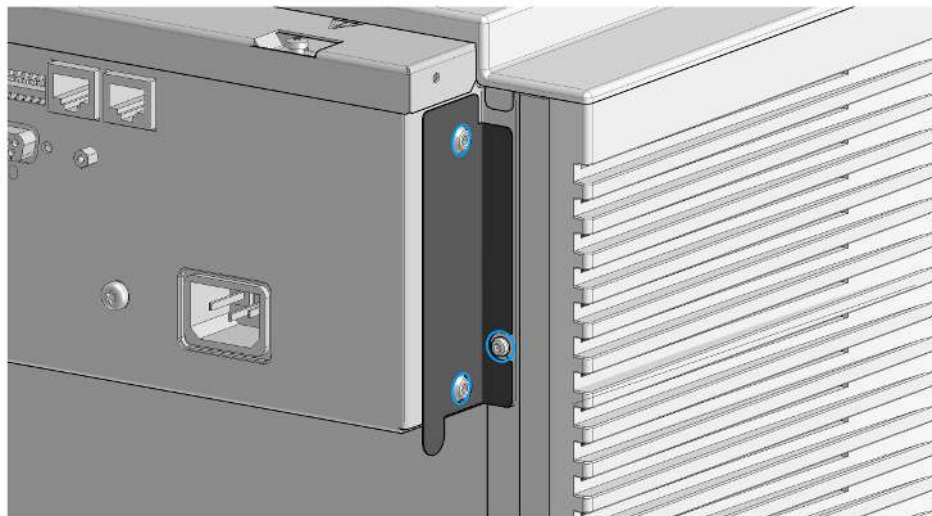


Install the Dust Filter Kit

- 9 Adhere the magnetic cover to the outer surface of the electronic box where the cooling fan is located. This area is on the right side when viewed from the back of the module.

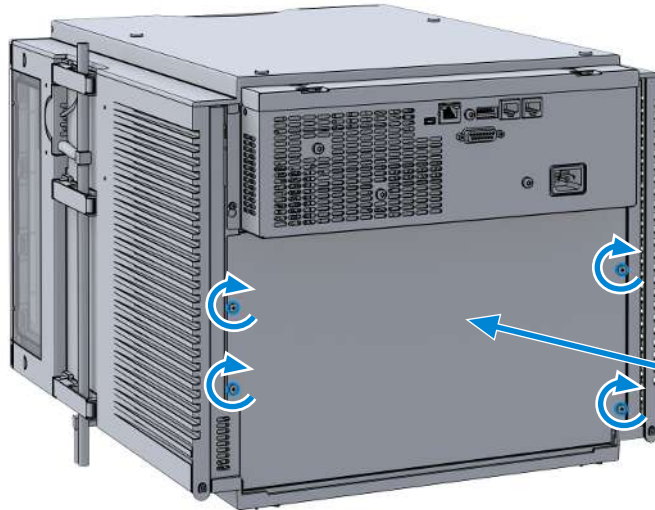


- 10 Ensure that the magnetic cover is oriented correctly and fully covers all venting holes.



11 Reinstall the sheet metal back cover and secure it with the four screws.

NOTE: If desired, reinstall the Sample Thermostat by following the procedure outlined in the module-specific user manual.



12 Reconnect the power cable to the module.

