

Agilent 1290 Infinity II Preparative Open-Bed Sampler/Collector Bottom Fume Hood kit

Technical Note

This technical note describes the Preparative Open-Bed Sampler/Collector Bottom Fume Hood kit.

Install the Bottom Fume Hood

The door stop of the bottom fume hood can be mounted on the left or right side of the Open-Bed Fraction Collector.

Parts required	p/n	Description
	G9321-68300	Bottom Fume Hood Kit

Preparations	
	<ul style="list-style-type: none">• Preparative Open-Bed Fraction Collector <i>connected to ventilation system</i>.• Parts unpacked• Protective film removed

WARNING

Improper installation

Malfunctioning or hazardous conditions

- Read the following instructions carefully.
- Only trained personnel is allowed to install the module to the lab venting systems.
- Respect the requirements as specified in this document.

WARNING

Insufficient ventilation

Organic solvent vapours can reach explosive concentrations when they are not removed sufficiently with an appropriate venting system.

- Connect the exhaust tube to the laboratory's venting system.
- Ensure a minimum air flow of 6 L/s through the exhaust tube.
- Ensure an underpressure in the venting system.



WARNING

Flammable solvent vapors

Fire or explosion hazard through electrostatic discharge

- Ground the venting tube, the analytical equipment and the venting system.
-

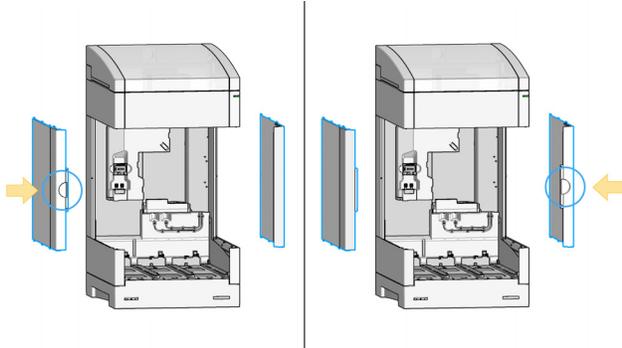
WARNING

Toxic, flammable and hazardous solvents, samples and reagents

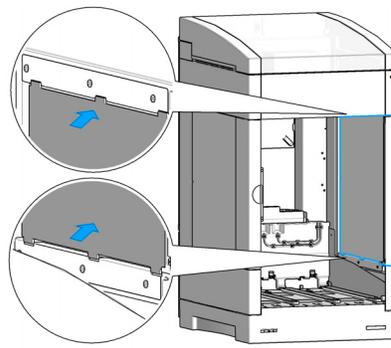
The handling of solvents, samples and reagents can hold health and safety risks.

- When working with these substances observe 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.
 - Do not use solvents with an auto-ignition temperature below 200 °C (392 °F). Do not use solvents with a boiling point below 56 °C (133 °F).
 - Avoid high vapor concentrations. Always keep the temperature in the sample compartment at least 25 K below the boiling point of the solvent used.
 - Do not operate the instrument in an explosive atmosphere.
 - Do not use solvents of ignition Class IIC according IEC 60079-20-1 (for example, carbon disulfide).
 - Reduce the volume of substances to the minimum required for the analysis.
 - Never exceed the maximum permissible volume of solvents (8 L) in the solvent cabinet. Do not use bottles that exceed the maximum permissible volume as specified in the usage guideline for solvent cabinet.
 - Ground the waste container.
 - Regularly check the filling level of the waste container. The residual free volume in the waste container must be large enough to collect the waste liquid.
 - To achieve maximal safety, regularly check the tubing for correct installation.
-

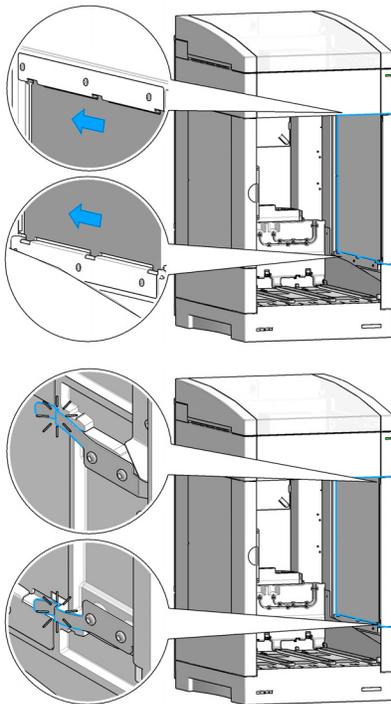
1 Decide on orientation of door opening (depends on installation of grip recess).



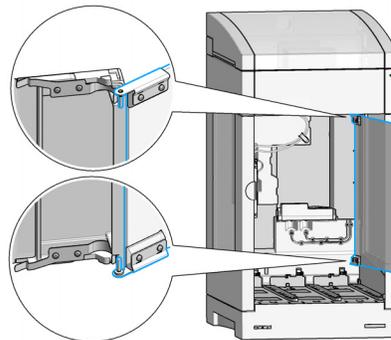
2 Place the brackets of the Bottom Fume Hood side panel into the gaps of the Fraction Collector side panel (bottom and top).



3 Push the side panel backwards until it clicks into place.

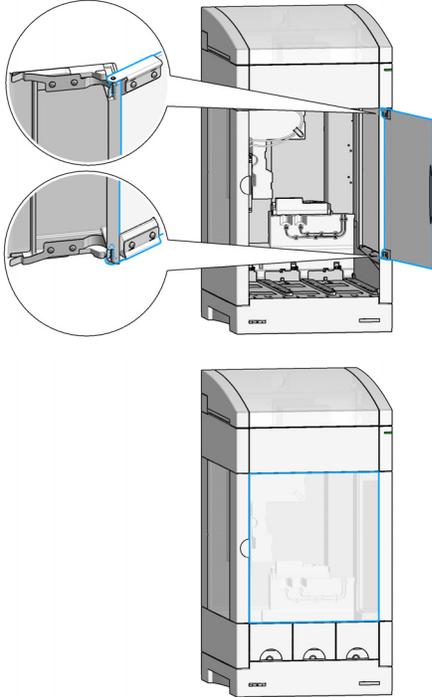


4 Position the front door with the two metal pins in direction to the side panel without grip recess.

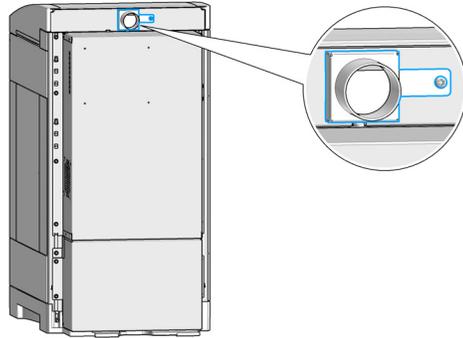


Install the Bottom Fume Hood

- 5** Press the metal pin into the spring clip to mount the door into the side panel (bottom and top).



- 6** Connect an exhaust hose (40 mm i.d.) to the top fume hood exhaust adapter and the ventilation system.



Next Steps:

- 7** Obtain and attach an exhaust tube (tube ID must be slightly greater than 40 mm) to the adapter. An exhaust tube is NOT provided with the Exhaust Tube Adapter kit. The exhaust tube must be mechanically stable, leak tight and resistant to the solvents used in your system.
- 8** Fix the exhaust tube to the adapter with an appropriate connector (e.g. hose clamp). Ground the exhaust tube.
- 9** The other outlet of the exhaust tube must be installed in a fume hood or connected to a vent system in a way, that the resulting air flow through the exhaust tube is at least 10 m³/h.



G7158-90200 Rev. B
SD-29000164
Edition: 10/2019

Printed in Germany

© Agilent Technologies, Inc 2018-2019

Agilent Technologies, Inc
Hewlett-Packard-Strasse 8
76337 Waldbronn
Germany