

Installation of the Infinity II Cooler/Thermostat Condensate Drainage Tubing Kit

Technical Note

This Technical Note describes the use and installation of the Infinity II Cooler/Thermostat Condensate Drainage Tubing Kit.

Delivery Checklist

Condensate Drainage Kit (5067-6208) contains:

Item #	Description
1	Condensate Collector
2	Single Drain Connector
3	Washer, ST, ID/OD 3.2/9 mm
4	Tubing, silicon rubber 2.16 m, ID/OD 6/9 mm (pre-cut)
5	Tubing connector, Y-shaped, ID 6.4 mm (with a 60 and 100 mm premounted tubing piece)
6	Tubing connector, 90°, ID 6.4 mm
7	Tubing holder clip, 19 x 19 x 8 mm

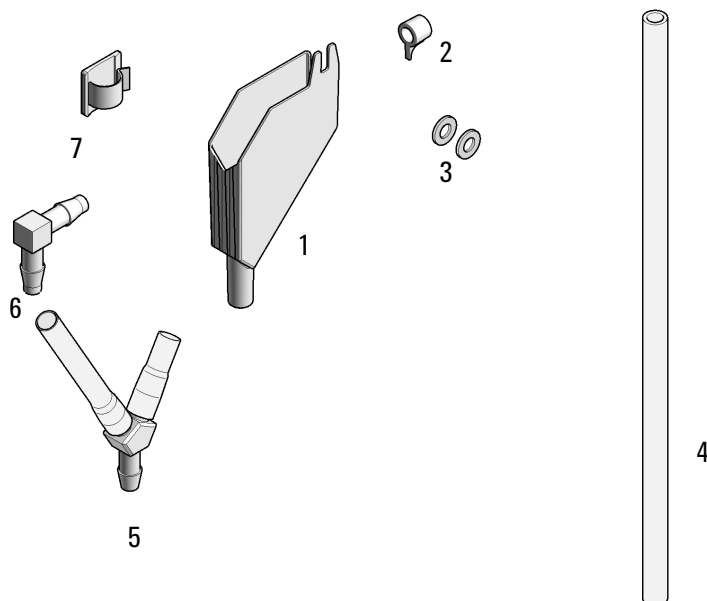


Figure 1 Content of Condensate Drainage Kit (5067-6208)

Installation of the Condensate Drainage Kit

Setup With the Condensate Collector Funnel for Bench Installations

Tools required	p/n	Description
	8710-1622	Torx screwdriver T15
OR	5023-3089	Torx key set

Parts required	#	Description
	1	Condensate Collector
	1	Single Drain Connector
	1	Washer, ST, ID/OD 3.2/9 mm
	1	Tubing, silicon rubber 2.16 m, ID/OD 6/9 mm (pre-cut)
	1	Tubing connector, 90 °, ID 6.4 mm

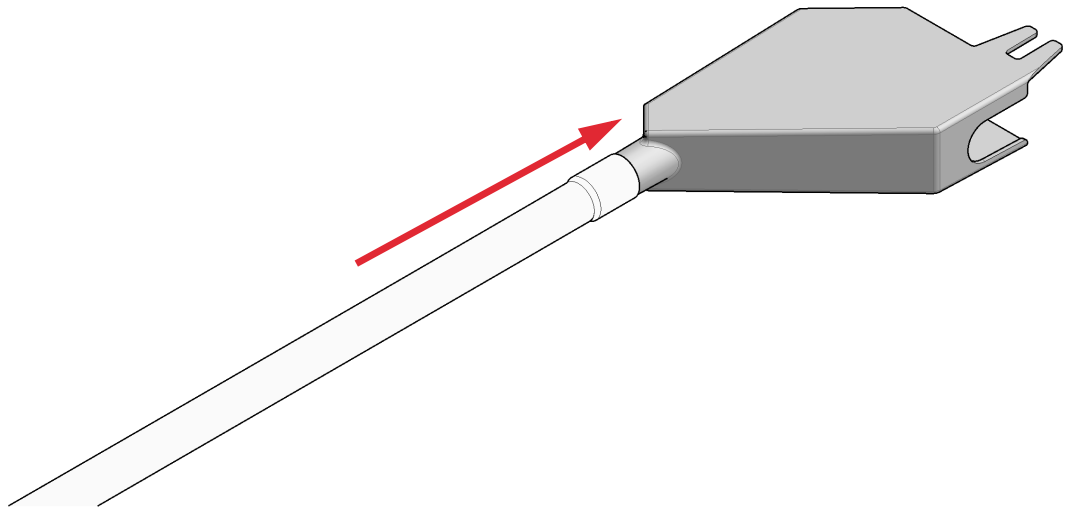
NOTE

The setup with the condensate collector funnel is suitable for bench installations only.

NOTE

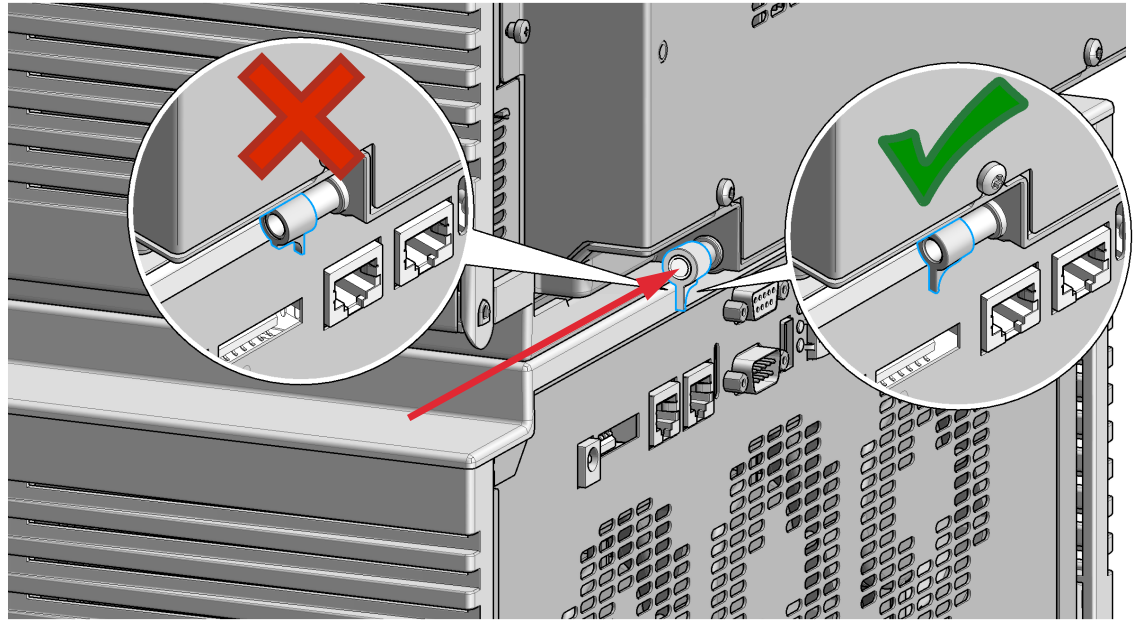
The condensate collector funnel is not suitable for Flex Bench systems or when the sampler is the lowermost module in a bench installation.

- 1 Attach the condensate tube to the outlet port of the condensate collector funnel.

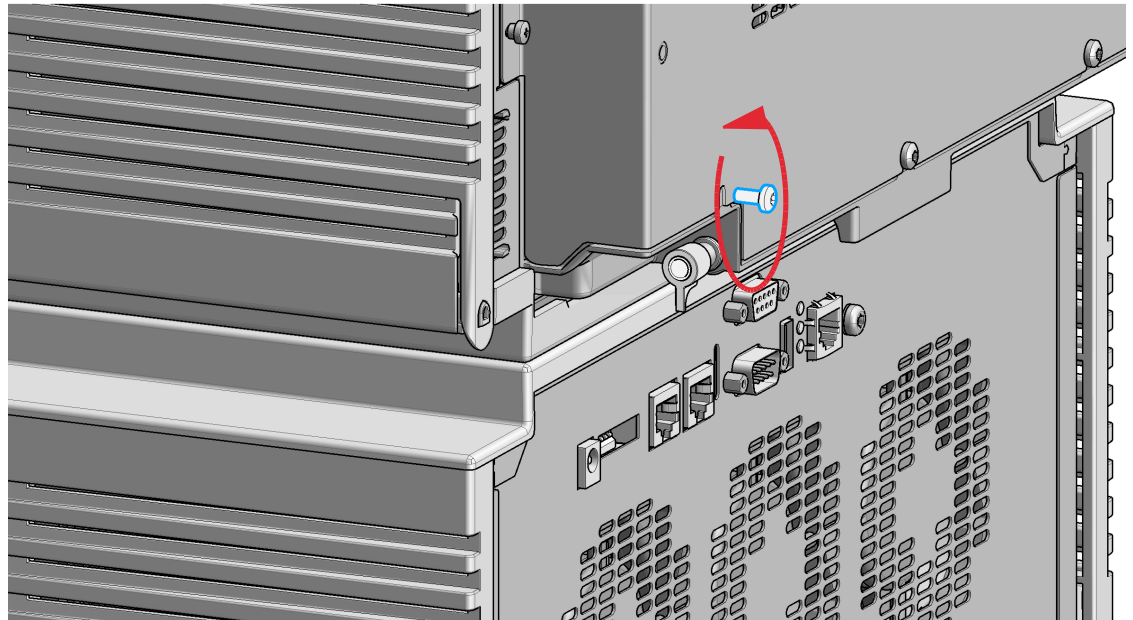


Installation of the Condensate Drainage Kit

- 2 Mount the drain connector on the condensate drainage outlet tube. Ensure the correct orientation of the spout.

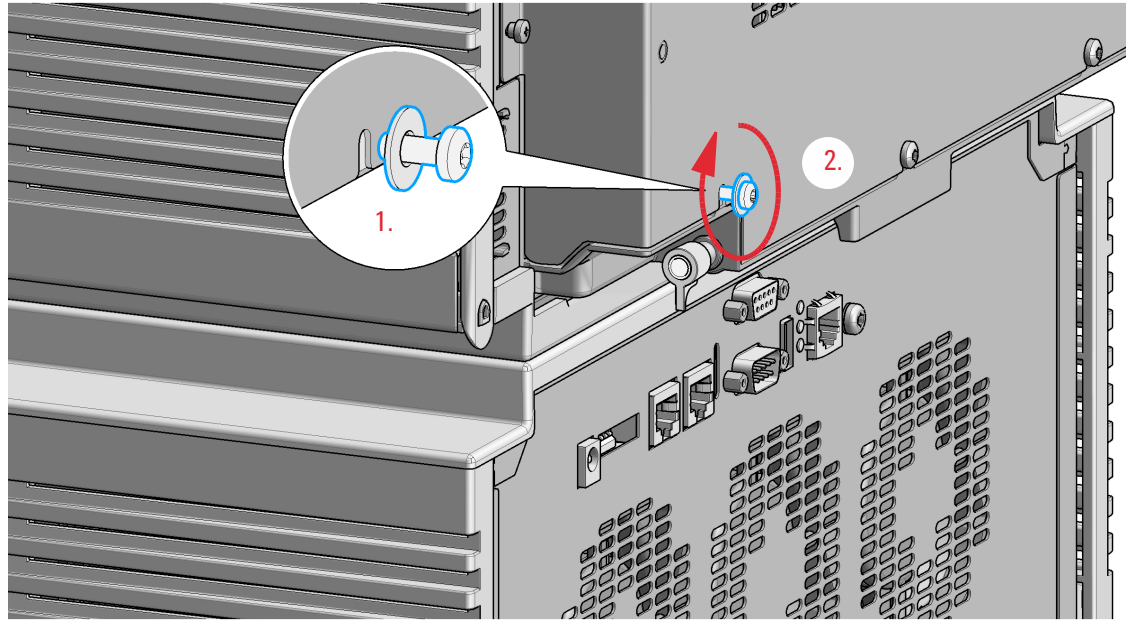


- 3 Remove the screw situated above the condensate drainage outlet tube.

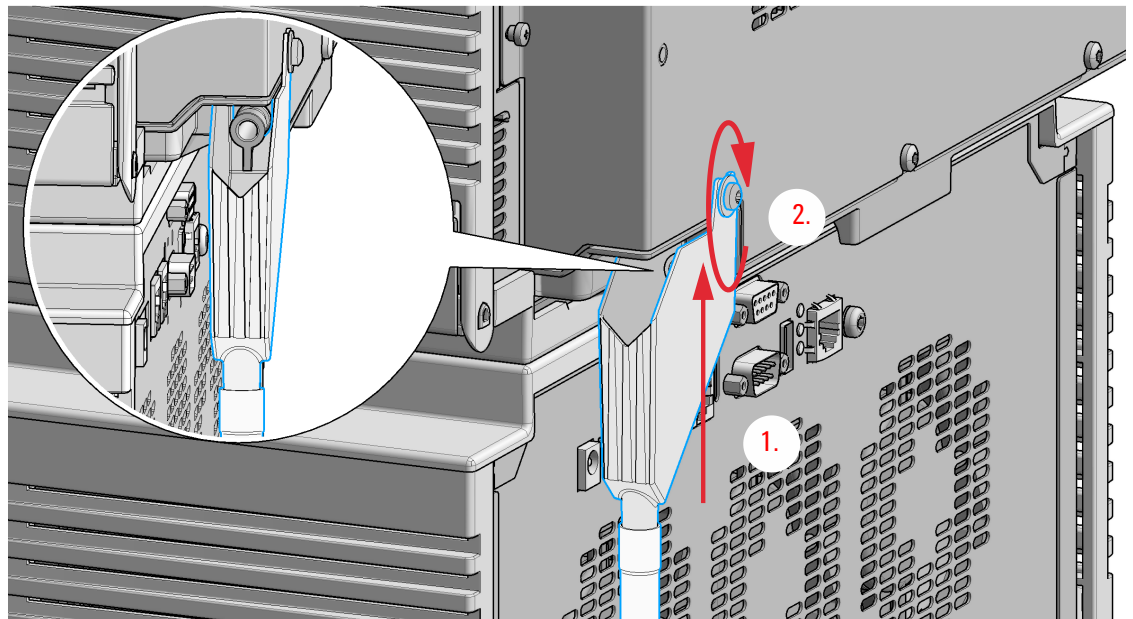


Installation of the Condensate Drainage Kit

- 4 Place the washer over the thread of the screw (1). Screw the screw and washer halfway into the hole in the back of the cooler/thermostat (2).

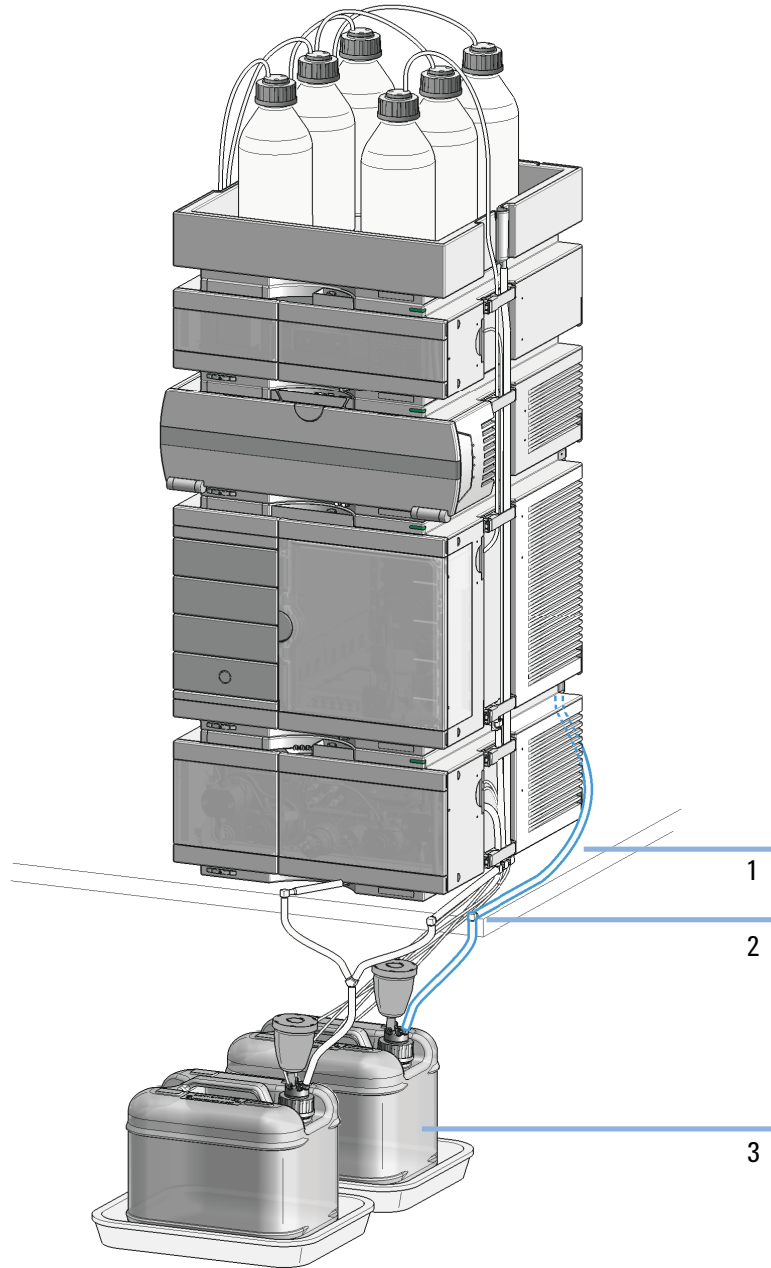


- 5 Position the condensate collector funnel underneath the condensate drainage outlet tube (1) and fix it to the back of the cooler/thermostat by tightening the screw (2). Ensure correct orientation and avoid over-tightening the screw.



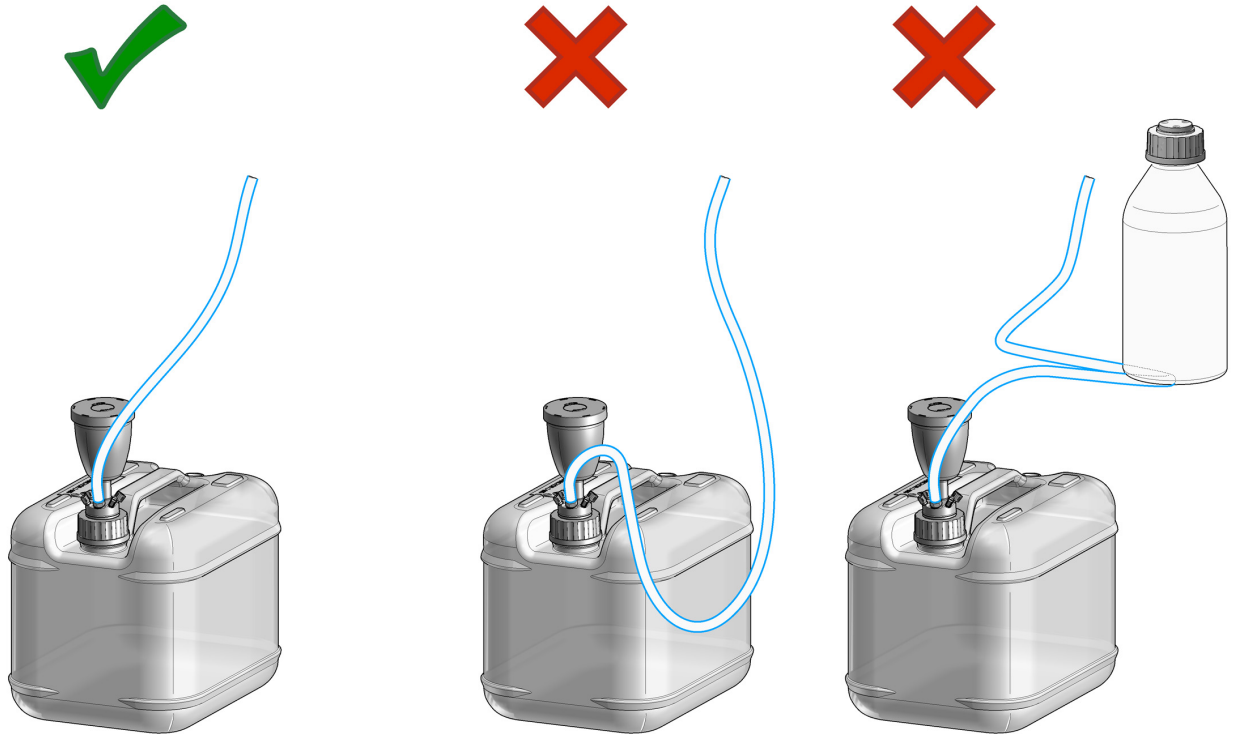
Installation of the Condensate Drainage Kit

- 6 Shorten the condensate tube so that it runs straight into the waste container without any unnecessary detour (1). If needed, use the 90 ° tubing connector provided in the kit to eliminate uphill sections, which might occur at the edge of the bench (2). Agilent recommends the use of a separate canister for condensate collection to avoid drainage problems (3).



Installation of the Condensate Drainage Kit

- 7 Ensure that the tubing runs straight into the waste canister without any bends or joints and it is not hindered by any mechanical obstacle. Agilent recommends using a 6 L waste canister equipped with a suitable InfinityLab Stay Safe cap for optimal condensate handling. If you decide to use your own waste solution, make sure that the tubes don't immerse in the liquid.



NOTE

Depending on the ambient conditions in the lab, the amount of condensate can vary from 200 mL to 2 L per day. Do not fill the waste container for the condensate to the top. Regularly empty the waste container.

Setup with the Y-piece Tubing Connector for Bench Installations and Flex Bench Systems

Parts required	Description
	Tubing connector, Y-shaped, ID 6.4 mm (with a 60 and 100 mm premounted tubing piece)
	Tubing holder clip, 19 x 19 x 8 mm
	Tubing, silicon rubber 2.16 m, ID/OD 6/9 mm (pre-cut)
	Tubing connector, 90 °, ID 6.4 mm

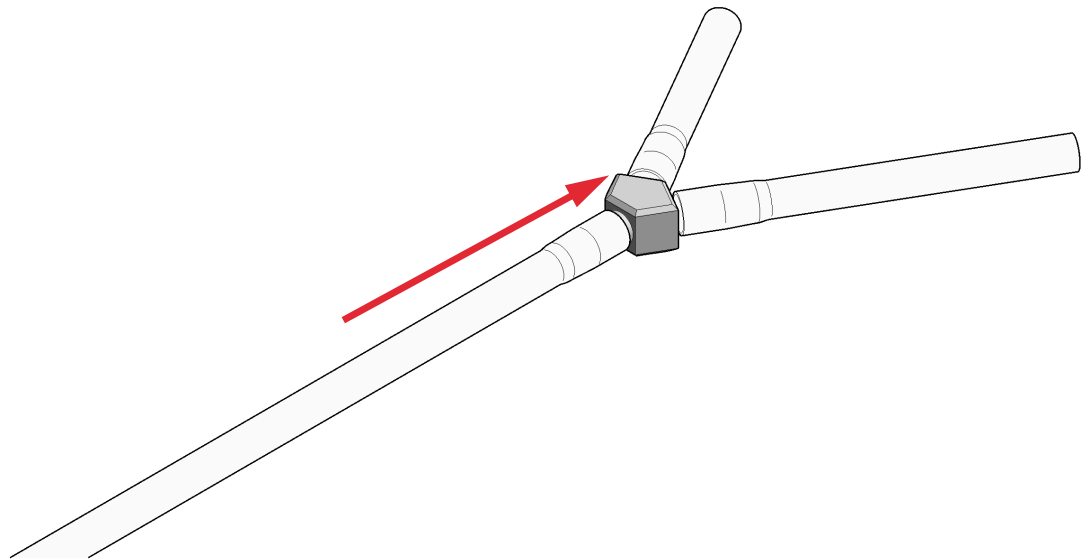
NOTE

The setup with the Y-piece tubing connector is equally suitable for bench installations and Flex Bench systems.

NOTE

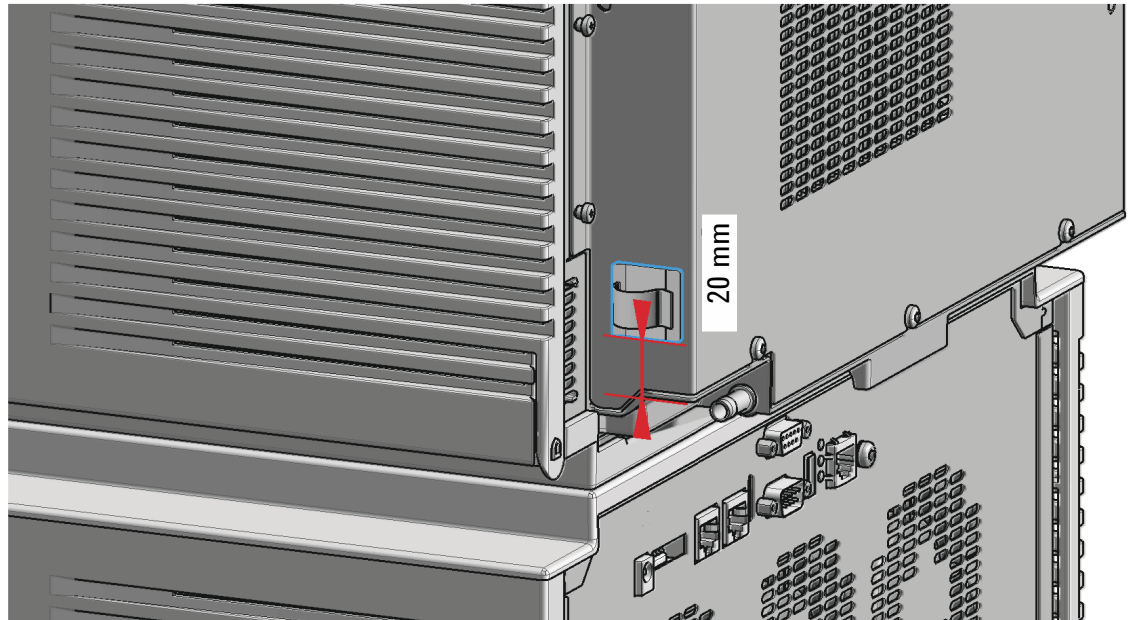
Agilent Infinity II autosamplers with a cooler/thermostat installed should not be used as the lowermost module in an HPLC stack to ensure adequate condensate drainage.

- 1 Attach the condensate tube to the free end of the Y-piece tubing connector.

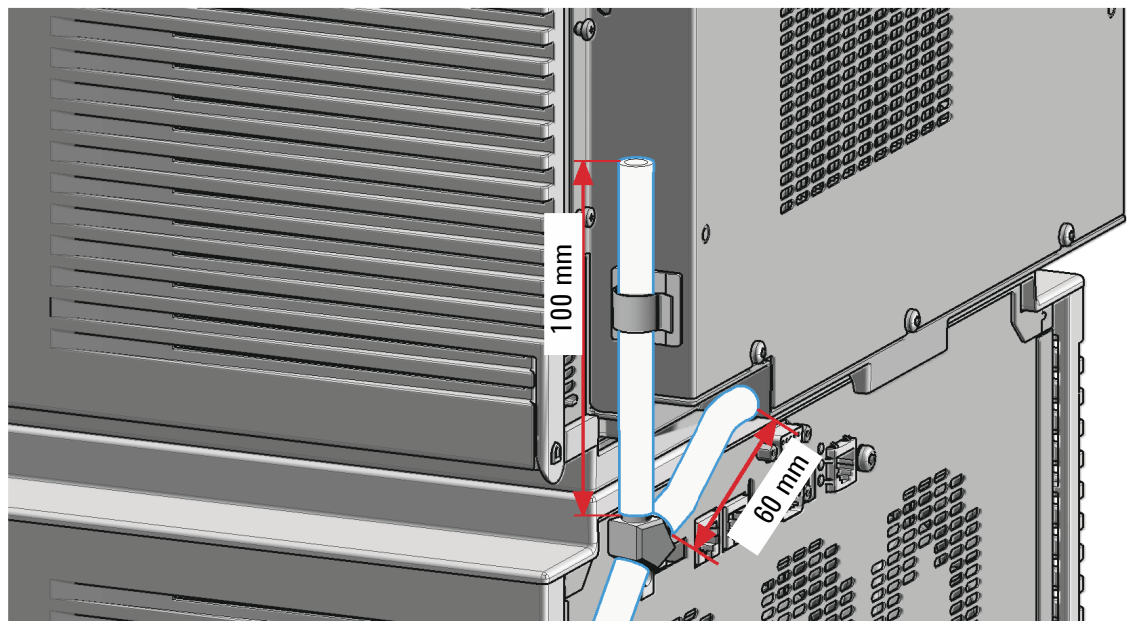


Installation of the Condensate Drainage Kit

- 2 Adhere a tubing holder clip to the side of the cooler/thermostat where the condensate drainage outlet tube is situated. Ensure a distance of 20 mm from the bottom edge.

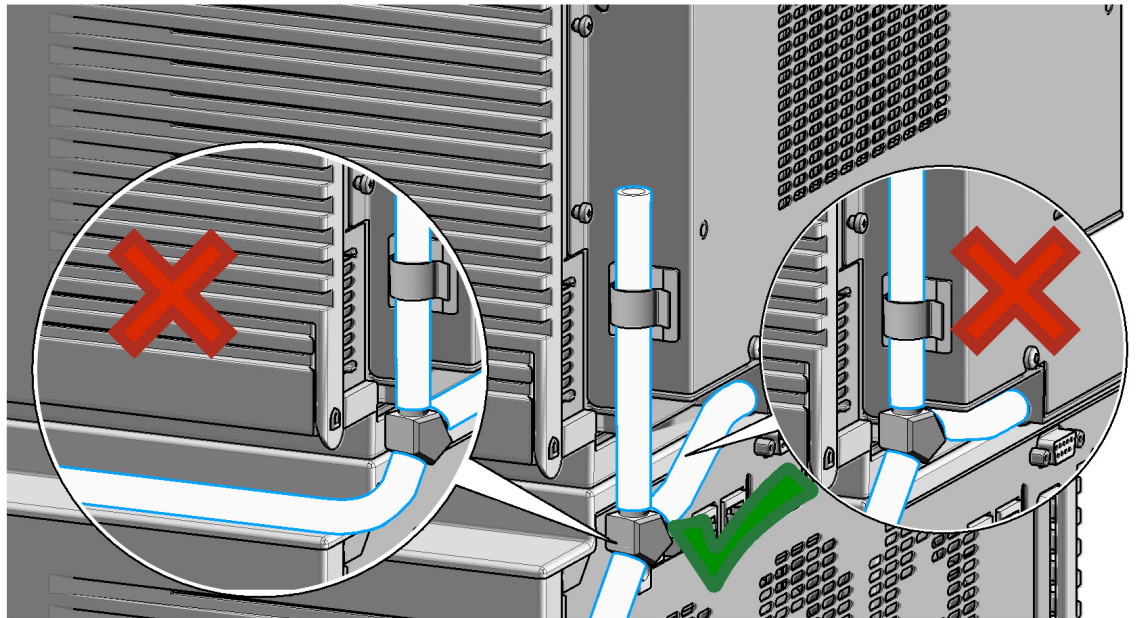


- 3 Mount the shortest tube (60 mm) of the condensate tubing with the Y-piece on the condensate drainage outlet tube and fix the venting tube (100 mm) in the tubing holder clip.



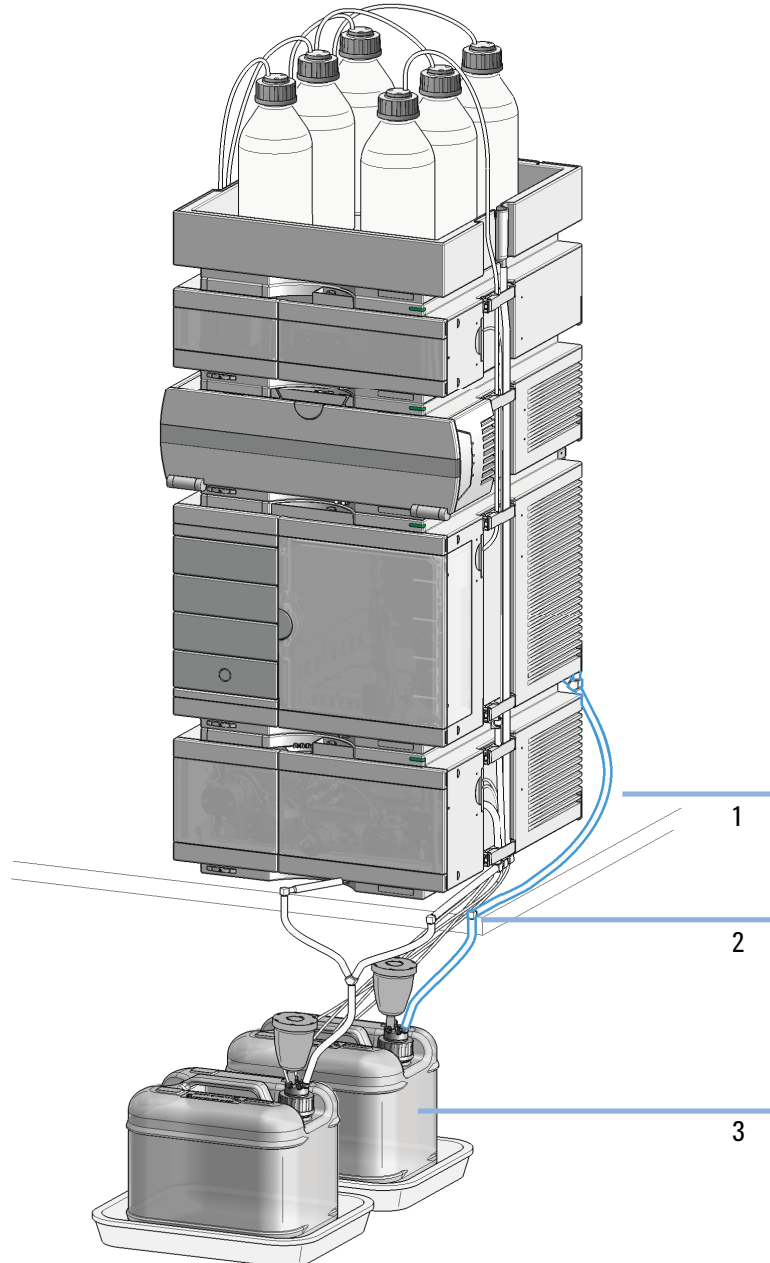
Installation of the Condensate Drainage Kit

- 4 Ensure that the condensate handling system is installed in a way that allows a continuous slope for the drained liquid. Horizontal or uphill sections may hinder the drainage.



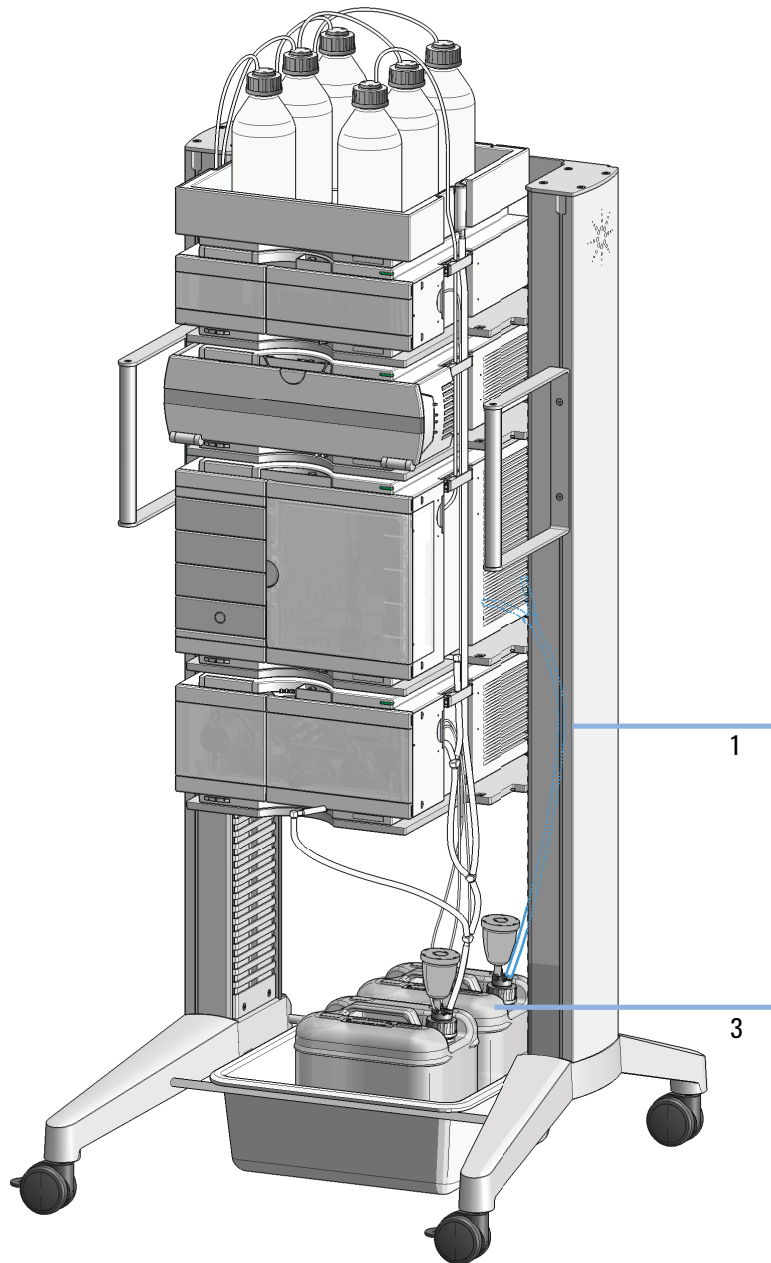
Installation of the Condensate Drainage Kit

- 5 Shorten the condensate tubing so that it runs straight into the waste container without any unnecessary detour (1). If needed, use the 90 ° tubing connector provided in the kit to eliminate up-hill sections, which might occur at the edge of the bench (2). Agilent recommends the use of a separate canister for condensate collection to avoid drainage problems (3).
 - a Routing the condensate tubing for bench installations.

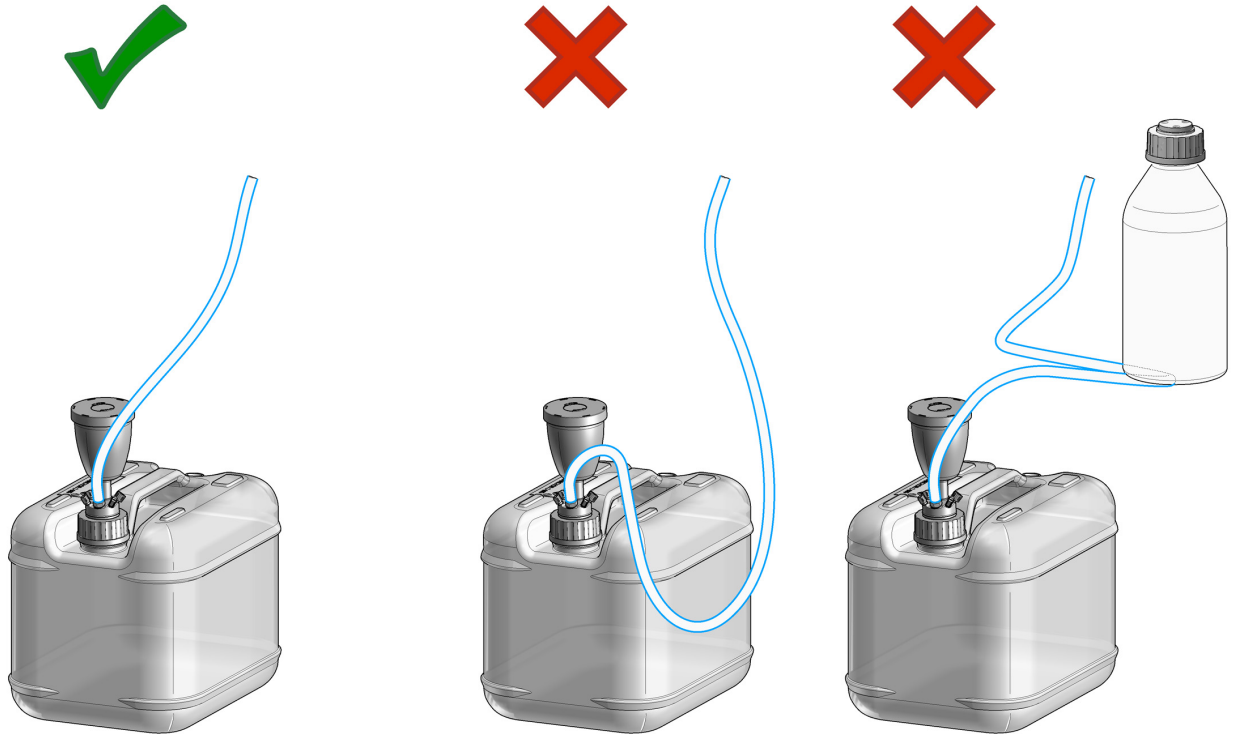


Installation of the Condensate Drainage Kit

- b Routing the condensate tubing for a Flex Bench system.



- 6 Ensure that the tubing runs straight into the waste canister without any bends or joints and it is not hindered by any mechanical obstacle. Agilent recommends using a 6 L waste canister equipped with a suitable InfinityLab Stay Safe cap for optimal condensate handling. If you decide to use your own waste solution, make sure that the tubes don't immerse in the liquid.



NOTE

Depending on the ambient conditions in the lab, the amount of condensate can vary from 200 mL to 2 L per day. Do not fill the waste container for the condensate to the top. Regularly empty the waste container.

