



# Agilent 1260 Infinity II Preparative/Analytical Fraction Collector Install Exhaust Tube Adapter Kit G1364-68709

## Technical Note

This technical note describes the installation of the Exhaust Tube Adapter Kit (G1364-68709) to one of the following modules:

- 1260 Infinity II Preparative Fraction Collector (G1364E),
- 1260 Infinity II Analytical Fraction Collector (G1364F),
- 1260 Infinity II Bio-inert Fraction Collector (G5664B).

## Exhaust Adapter Kit

The Exhaust Tube Adapter Kit (G1364-68709) contains:

p/n	Description
G1364-03260	Exhaust Tube Adapter
5067-5406	Screw Tx20, M4 x 8 mm

## Application and Safety Information

The exhaust tube adapter offers you the capability to connect your fraction collector to a laboratory's venting system, allowing the adaptation of a tube (not included) with an ID greater than 40 mm to your fraction collector.

### 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 10 m<sup>3</sup> / h through the exhaust tube.
- Ensure an underpressure in the venting system.



**NOTE**

An air flow through the vent tube greater than 10 m<sup>3</sup>/h reduces the cooling performance of a thermostatted instrument and increases the evaporation of solvents in your fraction vessels.

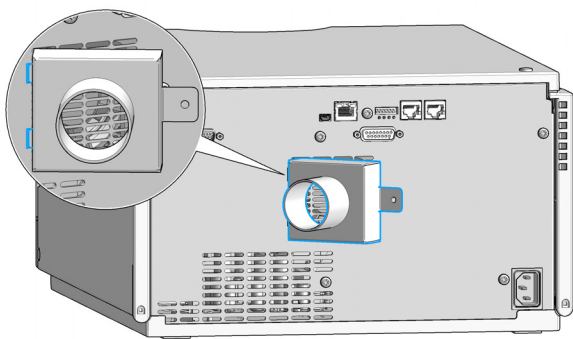
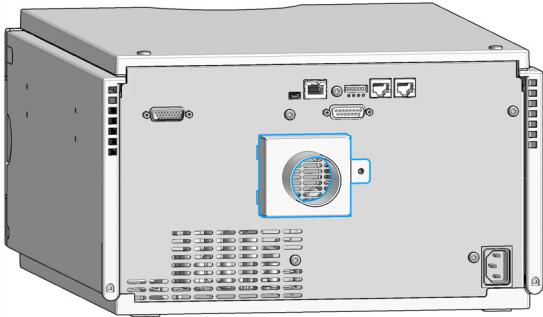
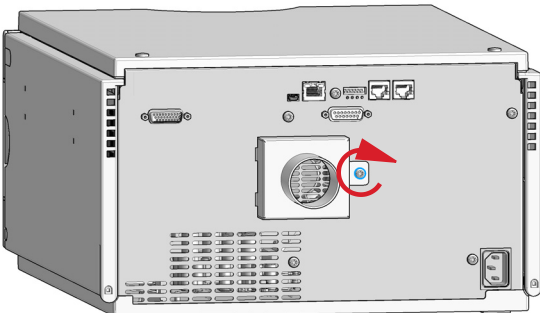
**WARNING**

**Flammable solvent vapors**

**Fire or explosion hazard through electrostatic discharge**

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

## Installation

<p><b>1</b> Attach the exhaust tube adapter to the left side slits of the rear fan exhaust grid with an angle of about 90 °.</p> 	<p><b>2</b> Push the adapter into the grid and turn it sideways, so that its latch covers the hole for the holding screw.</p> 
<p><b>3</b> Use the screw delivered in the Exhaust tube adapter kit to fasten the adapter latch to the rear of the fraction collector.</p> 	<p><b>Next Steps:</b></p> <p><b>4</b> 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.</p> <p><b>5</b> Fix the exhaust tube to the adapter with an appropriate connector (e.g. hose clamp). Ground the exhaust tube.</p> <p><b>6</b> 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 m<sup>3</sup>/h.</p>

