

www.agilent.com/chem/infinitylab-supplies

DE.3136689815

This information is subject to change without notice.

© Agilent Technologies, Inc. 2017, 2020

Published in US, March 5, 2020

Part Number 01200-90112

Agilent InfinityLab

Venting Valves

Engineered for efficiency and ease of use

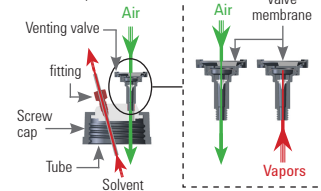


Venting Valve Information

The Agilent InfinityLab venting valve (p/n 5043-1190) plays an important role in maintaining your workspace, keeping it safe and clean. The venting valve blocks harmful vapors coming from the solvent bottle and assures the required ventilation while solvent is being pulled to the LC instrument. A filter membrane installed at the top of the valve avoids solvent contamination through dust particles. For optimum efficiency, a time strip installed on top of the filter indicates by when the filter has to be changed.

Venting Valve Working Principle

Membrane holds back dust and contaminant particles



Installation Procedure

1. Identify the connecting port on the cap. The venting valve can be installed on one of the connection ports below:



2. Install the venting valve on the connecting port and turn clockwise until you feel resistance.

3. Activate the time strip by pressing the bar with your finger.

The Time Strip

How to activate the time strip

Installed on top of the valve or the filter, a simple press with the finger on the bar starts the clock for a time of 6 months.



When Replacing the Filter or Valve

We recommend changing the InfinityLab venting valve (p/n 5043-1190) or the InfinityLab charcoal filter (p/n 5043-1193) when the red bar reaches the 6 months mark under the following typical conditions:

- 1 mL/min
- 5 days per week
- 8 hours per day
- Typical lab conditions at 20 °C

Note: The time strip is a visual reminder. It cannot provide real online information regarding the saturation and contamination of the filter and valve. The progress of the red bar is not linear over the 6 months period of time.