

Agilent InfinityLab LC Series

1260 Infinity II Delay Coil Organizer

Technical Note

This Technical Note describes the installation of the G9324A Delay Coil Organizer to an Agilent Infinity II system.

Contents

Delay Coil Organizer 2

Overview 2

Delivery Checklist 3

Parts 4

Installation 5

Assemble the Delay Coil Organizer 5

Assemble and Install the Leak Tray and Leak Sensor 7

Install the Delay Coil Organizer to another Module 8

Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column
Compartment 10



Agilent Technologies

Delay Coil Organizer

Overview

The Delay Coil Organizer allows the storage of up to three Delay Coil combinations to achieve the required delay volume, and to easily detect and dispose of any possible leaks.

The Delay Coil Organizer comes with a Pressure Relief Valve (PRV, as part of the accessory kit) for protection of the Fraction Collector inlet line.

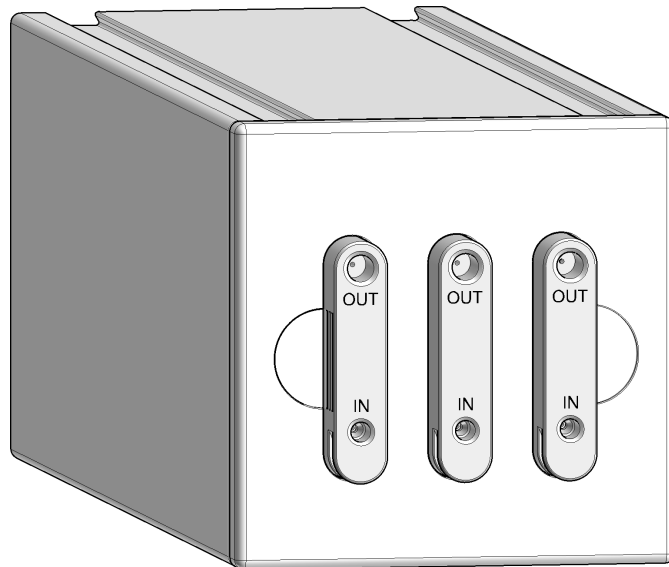


Figure 1 Delay Coil Organizer

Delivery Checklist

p/n	Description
G7163-60010	Delay Coil Box Assembly
G7163-68750	Accessory Kit Delay Coil Organizer
5067-6181	Knitted Delay Kit 1/2/5 m x 1.0 mm ESD Black (OPTIONAL)
5067-6800	Delay Coil Kit 4 – 8 mL/min ESD (OPTIONAL)
5067-6801	Delay Coil Kit 15 – 40 mL/min ESD (OPTIONAL)
5067-6802	Delay Coil Kit 40 – 80 mL/min ESD ¹ (OPTIONAL)
5067-6804	Delay Coil Kit 80 – 200 mL/min ESD ¹ (OPTIONAL)
5188-8053	ERI Remote / Leak Module (OPTIONAL)

¹ requires Adapter Nut FB 1/4-28-M 5/16-24-F ST (5023-3093), which comes as part of the Delay Coil Kit

NOTE

Delay coils are specified to 25 bar.

Accessory Kit

p/n	Description
5063-6527	Tubing assembly, i.d. 6 mm, o.d. 9 mm, 1.2 m (to waste)
5067-6655	Leak Sensor Assembly 3 m
5067-6857	PREP Pressure Relief Valve Kit
5067-5685	Clamp Guide Kit-IF-II
5043-0270	Leak plane
5043-0271	Holder leak plane
5043-0275	Clamp guide

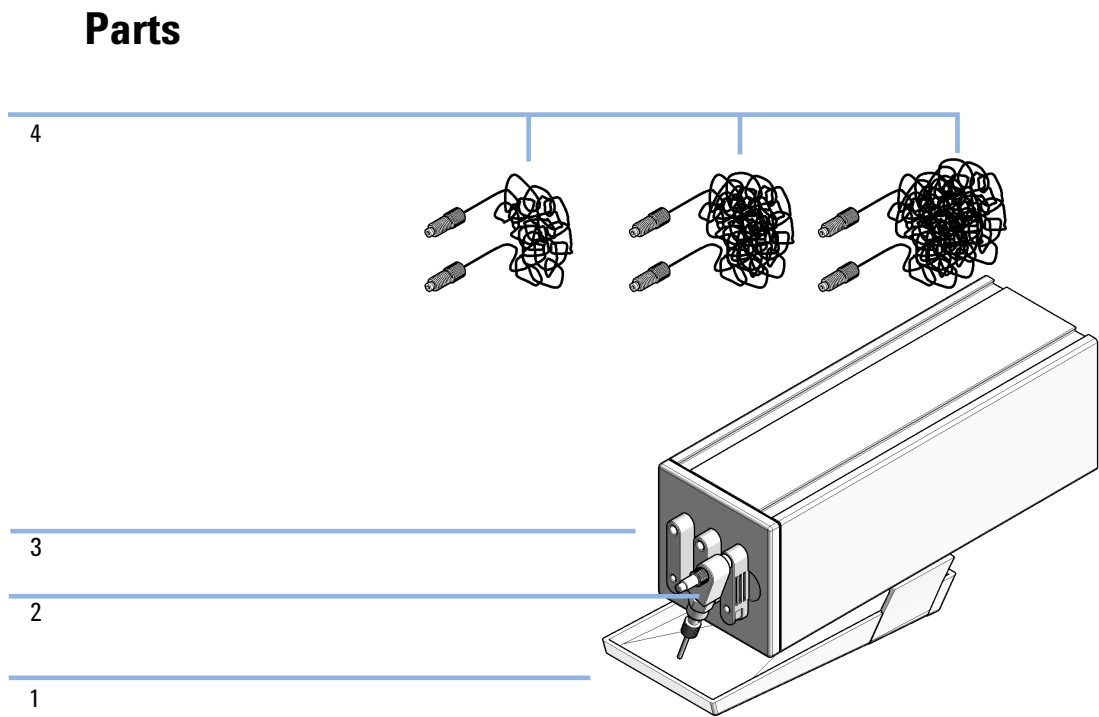


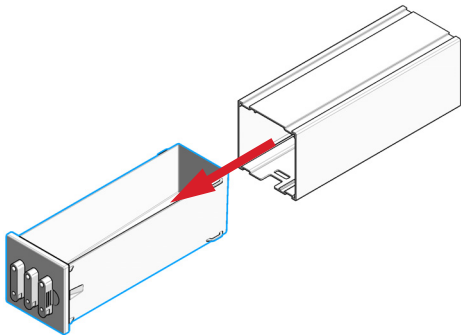
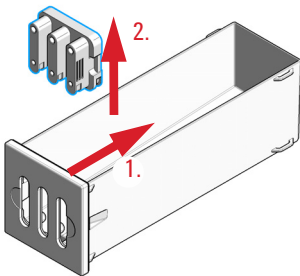
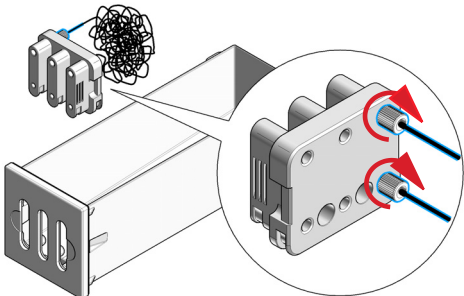
Figure 2 Delay Coil Organizer parts

Item	p/n	Description
1	5043-0270	Leak plane
2	5067-6857	PREP Pressure Relief Valve Kit
3	G7163-60010	Delay Coil Box Assembly
4	5067-6181	Knitted Delay Kit 1/2/5 m x 1.0 mm ESD Black (OPTIONAL)
	5067-6800	Delay Coil Kit 4 – 8 mL/min ESD (OPTIONAL)
	5067-6801	Delay Coil Kit 15 – 40 mL/min ESD (OPTIONAL)
	5067-6802	Delay Coil Kit 40 – 80 mL/min ESD (OPTIONAL)
	5067-6804	Delay Coil Kit 80 – 200 mL/min ESD (OPTIONAL)

Installation

You can install the Delay Coil Organizer either to the side of a particular module (see “[Install the Delay Coil Organizer to another Module](#)” on page 8), or in the 1290 Infinity II Preparative Column Compartment (see “[Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column Compartment](#)” on page 10).

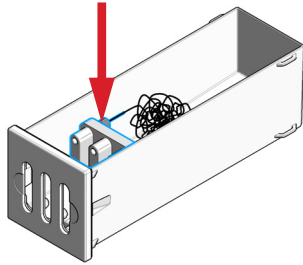
Assemble the Delay Coil Organizer

<p>1 Remove the delay coil organizer from its housing by pulling it to the front.</p> 	<p>2 Remove the distributor from the delay coil organizer.</p> 
<p>3 Install the Adapter Nut FB 1/4-28-M 5/16-24-F ST (5023-3093) for the Delay Coil Kit mL/min ESD (5067-6802) and the Delay Coil Kit mL/min ESD (5067-6804).</p> <p>4 Tighten the adapter nut with a torque of 2.3 Nm (about 10 ° after fingertight).</p>	<p>5 Mount the required delay coils to the available ports of the distributor, and fix them fingertight. Do not use any wrench or fitting handle.</p>  <div style="border: 1px solid black; padding: 5px; margin-top: 10px;"> <p>NOTE</p> <p>Use two vertical aligned ports for one coil (combination). You can install up to three combinations.</p> </div>

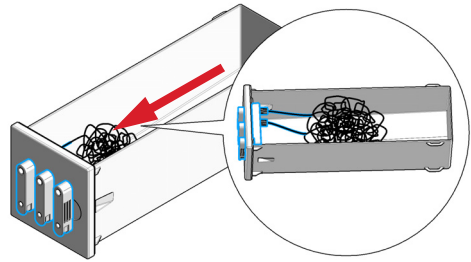
Installation

Assemble the Delay Coil Organizer

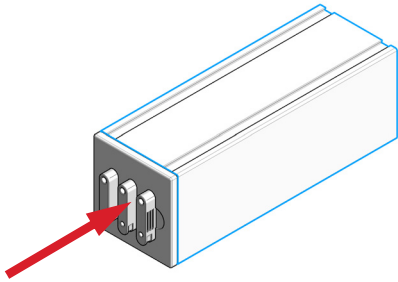
- 6** Place the distributor with the installed coil(s) back into the delay coil organizer.



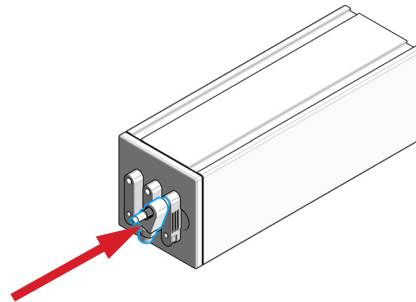
- 7** Install the distributor by pushing it through the holes in the module frame.



- 8** Slide the delay coil organizer into its housing to close it.

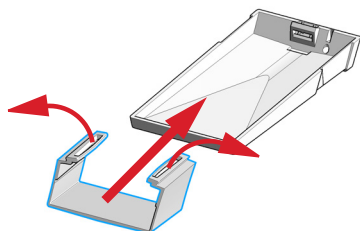


- 9** Install the Pressure Relief Valve (PRV) to the top port (OUT) of the Delay Coil (combination) to be used. This will indicate which port is in use.

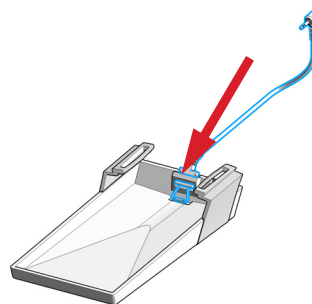


Assemble and Install the Leak Tray and Leak Sensor

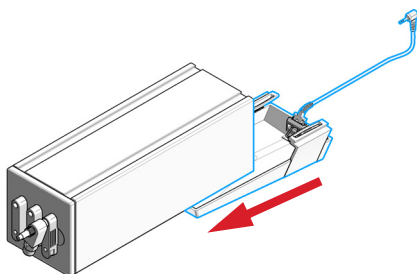
1 Attach the leak tray clamp to the leak tray.



2 Push the leak sensor into its holder.



3 Slide the leak tray onto the bottom of the Delay Coil Organizer.



Installation

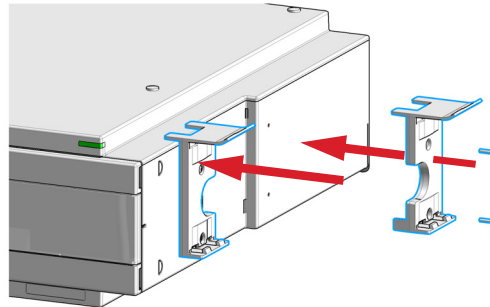
Install the Delay Coil Organizer to another Module

Install the Delay Coil Organizer to another Module

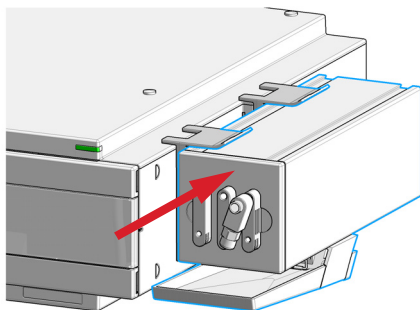
1 Assemble and install the leak tray and leak sensor to the delay coil organizer, see ["Assemble and Install the Leak Tray and Leak Sensor"](#) on page 7.

2 Ensure the power switch at the rear of the module is OFF and the power connector is unplugged.

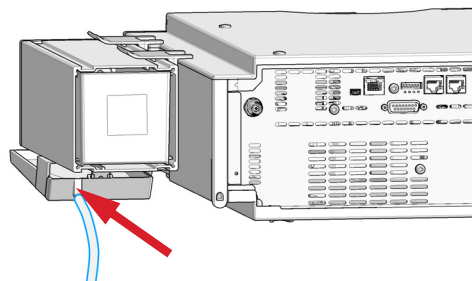
3 Identify the marks for the screws on the side panel of the cabinet assy and screw the clamp guides to the side panel of the cabinet.



4 Slide the delay coil organizer into the clamp guides from front to back.

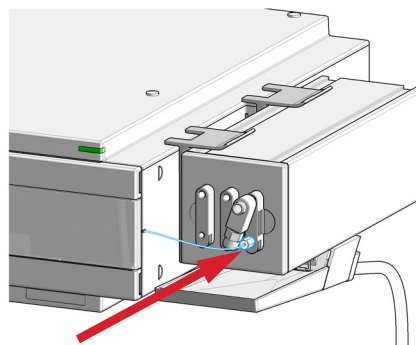


5 Attach the waste tubing to the leak plane and guide it to a proper waste container.

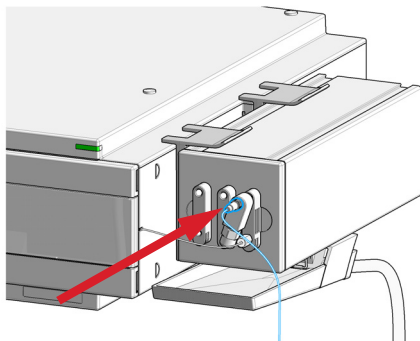


6 Connect the leak sensor cable to a module with ERI Remote connector (particularly G7159B, or G7161A/B), or to an ERI Remote / Leak Module (5188-8053).

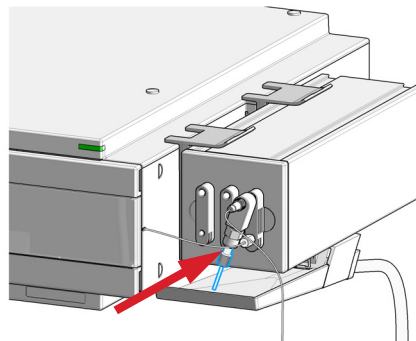
7 Install the capillary from the detector to the bottom port (IN) of the Delay Coil (combination) to be used.



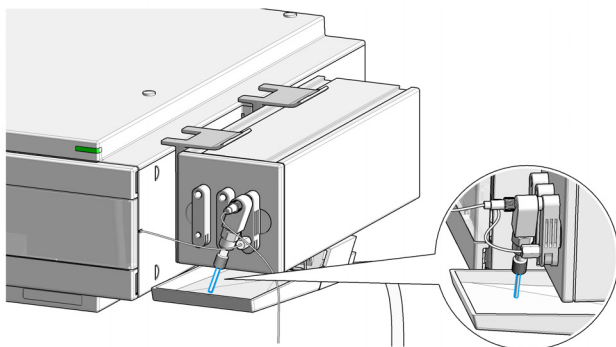
- 8** Install the tubing that goes to the fraction collector to the extended top front port of the PRV. Use the Fitting Handle to apply the correct torque.



- 9** Fix the waste tubing (1.6 mm x 650 mm ESD) to the bottom of the PRV.



- 10** Twist the PRV, such that the bottom does not get in the way of the bottom Delay Coil connection, but still ends inside the Leak Tray.



NOTE

All fittings used in combination with PRV, or inside the Delay Coil Organizer are $\frac{1}{4}$ -28 FB. Fittings to the bottom port of the Delay Coil Organizer are 10-32 coned.

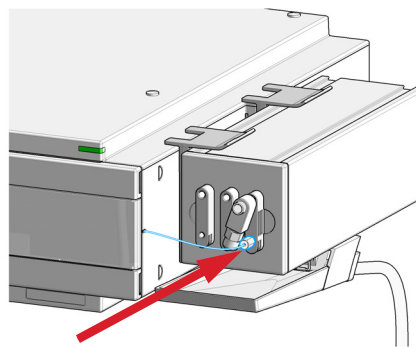
Installation

Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column Compartment

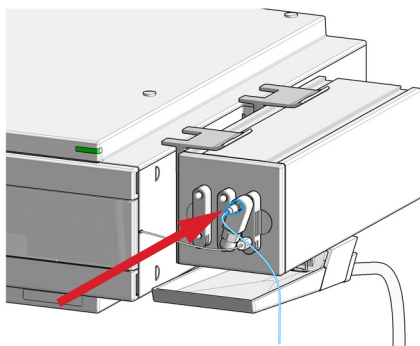
Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column Compartment

- 1** Install the Delay Coil Organizer into the top valve position of the 1290 Infinity II Preparative Column Compartment according to the Purification Capillary Kits Technical Note.

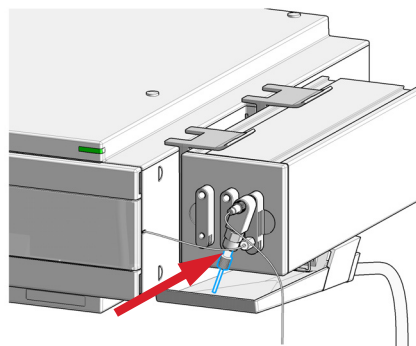
- 2** Install the capillary from the detector to the bottom port (IN) of the Delay Coil (combination) to be used.



- 3** Install the tubing that goes to the fraction collector to the extended top front port of the PRV. Use the Fitting Handle to apply the correct torque.

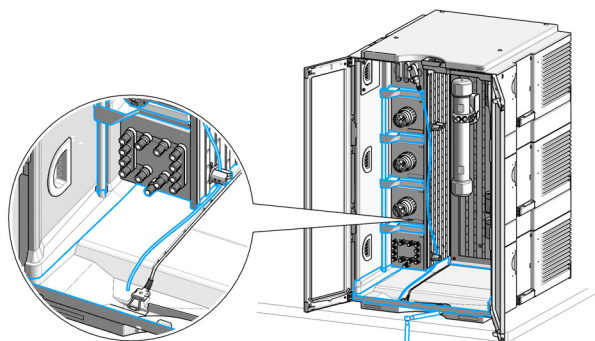


- 4** Fix the waste tubing (1.6 mm x 650 mm ESD) to the bottom of the PRV.



Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column Compartment

- 5** Twist the PRV, such that the bottom doesn't get in the way of the bottom Delay Coil connection, while the 650 mm waste line is guided alongside the small mounting plate inside the Column Compartment, to the Leak Plate and the tubing is fixed directly under the leak sensor assembly.

**NOTE**

All fittings used in combination with PRV, or inside the Delay Coil Organizer are $\frac{1}{4}$ -28 FB. Fittings to the bottom port of the Delay Coil Organizer are 10-32 coned.

Installation

Install the Delay Coil Organizer into the 1290 Infinity II Preparative Column Compartment



G9324-90000 Rev. B
SD-29000217 Rev.B
Edition: 10/2019

Printed in Germany

© Agilent Technologies, Inc 2017-2019

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