

Agilent InfinityLab LC Series  
1290 Infinity III LC System  
**System Manual**



# Notices

## Document Information

The information in this document also applies to 1260 Infinity II and 1290 Infinity II modules.

Document No: SD-29000160 Rev. E.01  
Edition: 05/2025

## Copyright

© Agilent Technologies, Inc.  
2014-2025

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

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

## Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

### WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

# Contents

## **In This Book 4**

### **1 Introduction 5**

Product Description of the 1290 Infinity III LC System 6

Features of the 1290 Infinity III LC System 7

System Components 8

Solutions 18

### **2 Installation 20**

Installing the System Modules 21

Optimizing the Stack Configuration 22

Integration Into the Network 30

Standard Capillary Setup 31

Handling Leak and Waste 33

### **3 Configuration Settings 45**

General Information on LAN Configuration 46

Instrument Configuration 47

Lab Advisor 49

### **4 Quick Start Guide 54**

Best Practices 55

Prepare a Run 59

Installation Checkout (Customer presence recommended) 69

### **5 Parts and Consumables 83**

HPLC System Tool Kit 84

InfinityLab Quick Connect and Quick Turn Fittings 85

Safety Caps and Solvent Bottles 90

InfinityLab Flex Bench Family 95



## In This Book

This manual covers the Agilent 1290 Infinity III LC System.

This chapter gives an introduction to the Agilent 1290 Infinity III LC System, the underlying concepts and features.

## **Product Description of the 1290 Infinity III LC System 6**

### **Features of the 1290 Infinity III LC System 7**

#### **System Components 8**

Product Description of the 1290 Infinity III Flexible Pump (G7104A) 9

Product Description of the 1290 Infinity III High-Speed Pump (G7120A) 10

Product Description of the 1290 Infinity III Multisampler (G7167B) 11

Product Description of the 1290 Infinity III Hybrid Multisampler (G7137B) 12

Product Description of the 1290 Infinity III Vialsampler (G7129B) 13

Product Description of the 1290 Infinity III Multicolumn Thermostat (G7116B) 14

Product Description of the 1290 Infinity III Diode Array Detector (G7117B) 15

Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B) 16

Product Description of the InfinityLab Level Sensing (G7175A) 17

#### **Solutions 18**

2D-LC Solution 18

High Dynamic Range DAD Solution 18

Method Transfer Solution 19

## Product Description of the 1290 Infinity III LC System

The Agilent 1290 Infinity III LC System has a high reliability and robustness, and improves analytical, instrument, and laboratory efficiency.

Its separation and detection performance delivers analysis data of high quality - the perfect front-end for high-end LC/MS systems.

Highest throughput for any application is achieved through high sample capacity and fast injection cycles, combined with new levels of usability.

Additionally, seamless integration in existing infrastructure and smooth method transfer from legacy equipment enable non-disruptive transition to highest productivity and lowest cost of ownership.

Designed for use with InfinityLab Assist, Level Sensing and Sample ID Reader.

## Features of the 1290 Infinity III LC System

- More chromatographic resolution: specially designed components in the sample flow path achieve low system dispersion.
- Higher peak capacity for challenging separations: switch easily between single dimension UHPLC and 2D-LC.
- Lower carryover for uncompromised data quality: multiwash capabilities of the Multisampler reduce carryover to less than 9 ppm even for challenging compounds such as chlorohexidine.
- Unique detection capabilities: combine low detection limits with an ultra-wide dynamic range by using the new 1290 Infinity III HDR-DAD or 1290 Infinity III ELSD.
- Faster injection cycles with dual needle injection for higher sample throughput.
- Higher sample capacity per bench space: up to 6144 samples within the footprint of a standard Agilent stack.
- Significantly better usability: dead-volume-free UHPLC fluidic connections can be easily achieved with the new InfinityLab Quick Connect fittings.
- Flexibility for all applications: due to the wide power, temperature and automatically scalable injection range, gradient options and intelligent system emulation technology.
- Seamless transfer of methods between LCs, regardless of the brand – facilitated by Intelligent System Emulation Technology (ISET) delivering unchanged retention time and peak resolution.
- Seamless integration in your chromatography data system: Agilent's Instrument Control Framework (ICF) enables smooth control of Agilent LC instrumentation through third-party chromatography data systems.
- Equipped with InfinityLab Assist - adds an Intuitive User Interface, Automated Workflows, Predictive Maintenance and Assisted Troubleshooting.

## System Components

The Agilent 1290 Infinity III LC System can be set up with two different types of pumps: with the G7104A Flexible Pump, combining the performance of a high-pressure mixing UHPLC pump with the flexibility of a low-pressure mixing UHPLC pump, or with the G7120A High-Speed Pump, allowing you to run fast gradients for high laboratory efficiency.

The Agilent 1290 Infinity III LC System consists of the following components:

- Flexible Pump (G7104A) or High-Speed Pump (G7120A)
- Multisampler (G7167B), Hybrid Multisampler (G7137B) or Vialsampler (G7129B)
- Multicolumn Thermostat (G7116B)
- Diode Array Detector (G7117B) or Variable Wavelength Detector (G7114B)
- Solvent Cabinet or InfinityLab Level Sensing (G7175A)

The Agilent 1290 Infinity III LC System is described in more detail in the following sections. All modules are stackable, see [Optimizing the Stack Configuration](#) on page 22.

For specifications, please refer to the individual module user documentation.



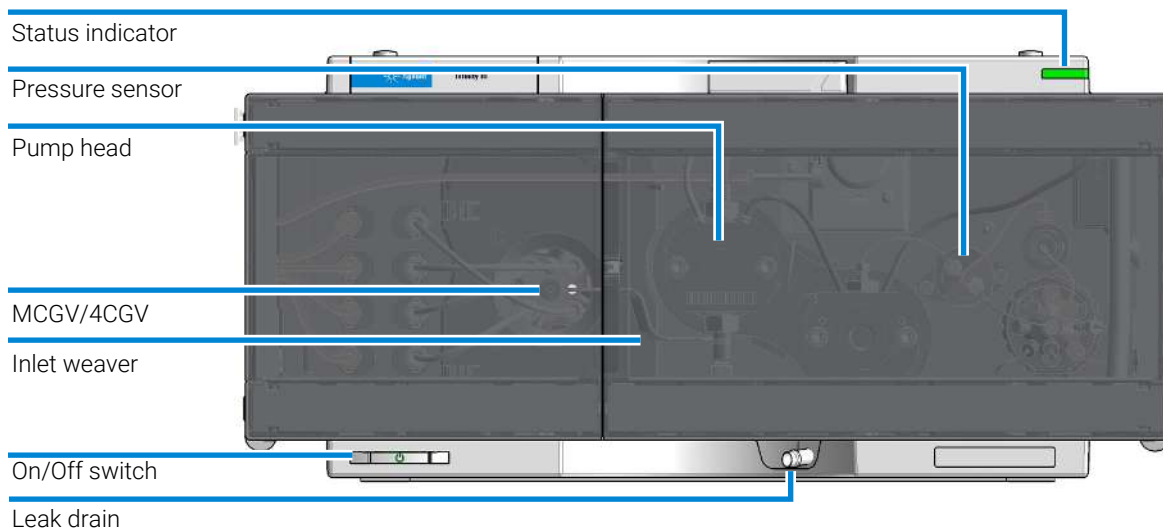
## Product Description of the 1290 Infinity III Flexible Pump (G7104A)

The Agilent 1290 Infinity III Flexible Pump improves your efficiency by combining the performance of a high-pressure mixing UHPLC pump with the flexibility of a low-pressure mixing UHPLC pump.

The new 1290 Infinity III LC power range has a high instrument efficiency, allowing you to run any HPLC and UHPLC method. ISET enables you to transfer existing methods from different instrument modules - current Agilent systems as well as instruments from other manufacturers.

The Agilent Inlet Weaver mixer, active damping or the optional Agilent Jet Weaver mixers for additional mixing capacity achieve high analytical efficiency.

The established multipurpose valve enhances laboratory efficiency by adding useful functionalities, for example, mixer in/out switch, filter backflush or automatic purge, and BlendAssist software simplifies your workflow with accurate buffer/additive blending.



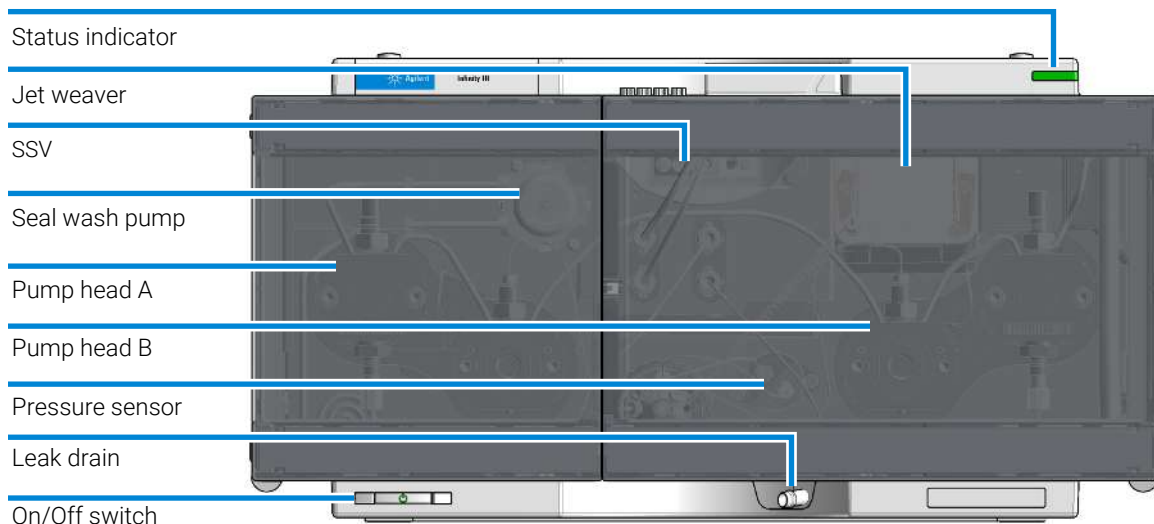
**Figure 1:** Overview of the Flexible Pump

## Product Description of the 1290 Infinity III High-Speed Pump (G7120A)

The Agilent 1290 Infinity III High-Speed Pump can enhance your efficiency through high speed and chromatographic performance. This binary pump is designed for full flexibility when running HPLC and UHPLC applications using narrow-bore columns and sub-2  $\mu\text{m}$  particles.

The 1290 Infinity III High-Speed Pump power range combines ultrahigh pressure up to 1300 bar and high analytical flow rates up to 5 mL/min for maximum UHPLC performance. Exceptionally low delay volumes down to 10  $\mu\text{L}$  and high flow rates enable fast gradients over the entire composition range from 1 to 99% for shortest run times and maximum resolution.

Highest retention time reproducibility and accuracy are achieved through independently controlled high-definition pump drives and active pressure damping.



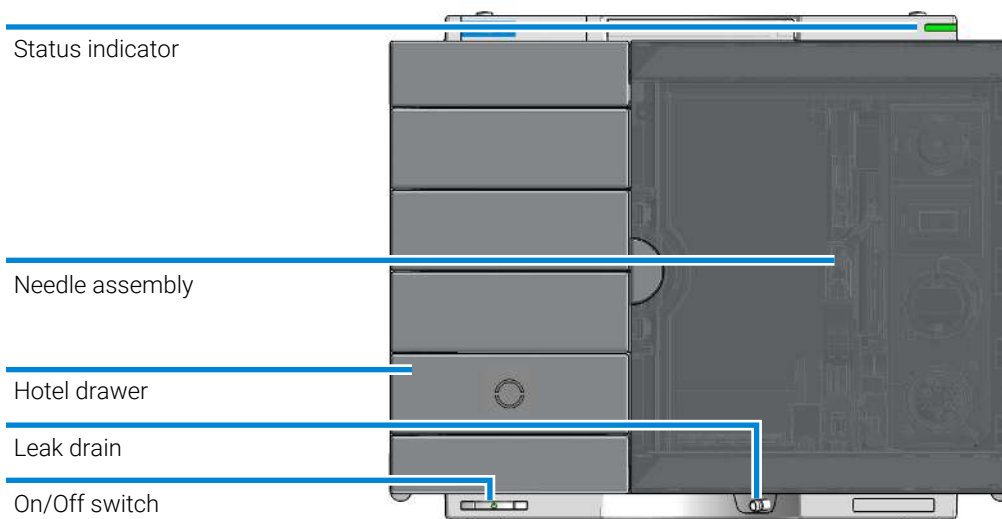
**Figure 2:** Overview of the High-Speed Pump

## Product Description of the 1290 Infinity III Multisampler (G7167B)

The Agilent 1290 Infinity III Multisampler can handle both vials and microtiter plates with ease and efficiency up to 1300 bar system pressure, optimized on chromatographic performance.

In fact, this compact module has the capacity to house up to 6144 samples, all inside the Agilent stack footprint and the robotics to smoothly inject each into the chromatograph in turn.

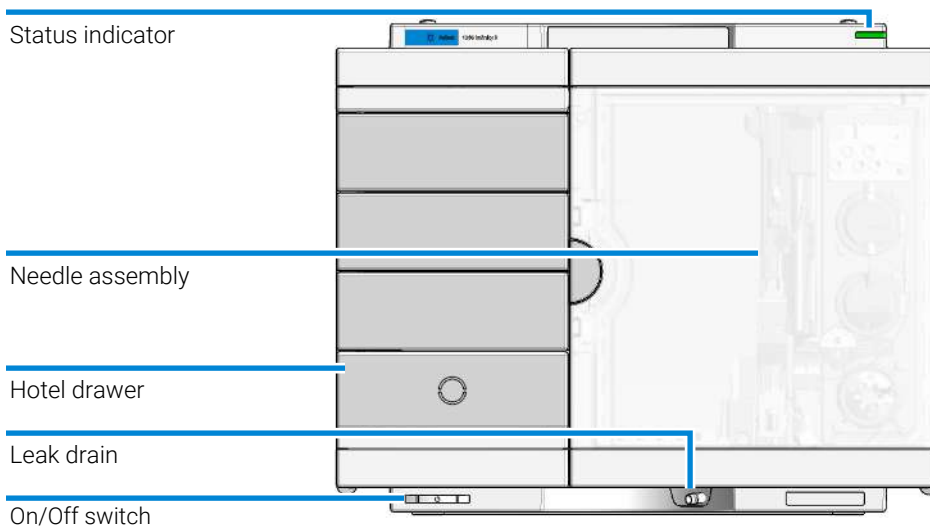
With the multi-wash capability, you can reduce carryover to less than 9 parts per million.



**Figure 3:** Overview of the Multisampler

## Product Description of the 1290 Infinity III Hybrid Multisampler (G7137B)

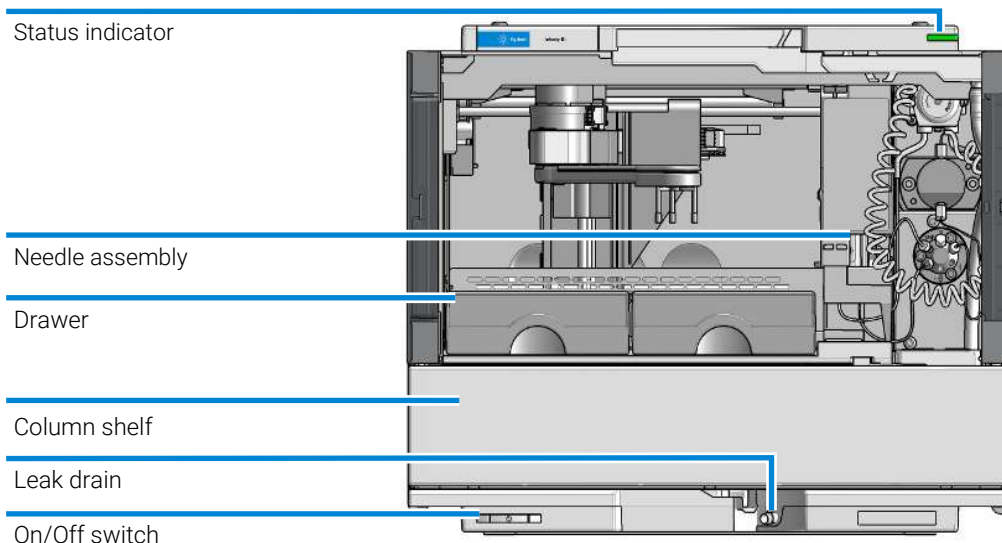
The Agilent 1290 Infinity III Hybrid Multisampler offers both standard flow-through and Agilent Feed Injection modes. For temperature-sensitive samples, a compressor-based thermostating device can be added. It can handle both vials and microplates with ease and efficiency up to 1300 bar system pressure, being optimized for high flexibility. This compact module can house up to 6144 samples, all inside the Agilent stack footprint and with the robotics to inject each into the chromatograph in turn. Its biocompatible sample flow path minimizes unwanted surface interaction of the sample.



**Figure 4:** Overview of the Hybrid Multisampler

## Product Description of the 1290 Infinity III Vialsampler (G7129B)

The Agilent 1290 Infinity III Vialsampler is an autosampler designed for UHPLC applications up to 1300 bar. It provides the reliability, safety, and ease-of-use needed for routine pharmaceutical tasks and quality control, as well as for environmental and food analyses. It can house optionally the integrated column compartment for two LC columns with temperature control up to 80 °C as well as a sample thermostat for stable temperatures from 4 °C to 40 °C, all within one module.



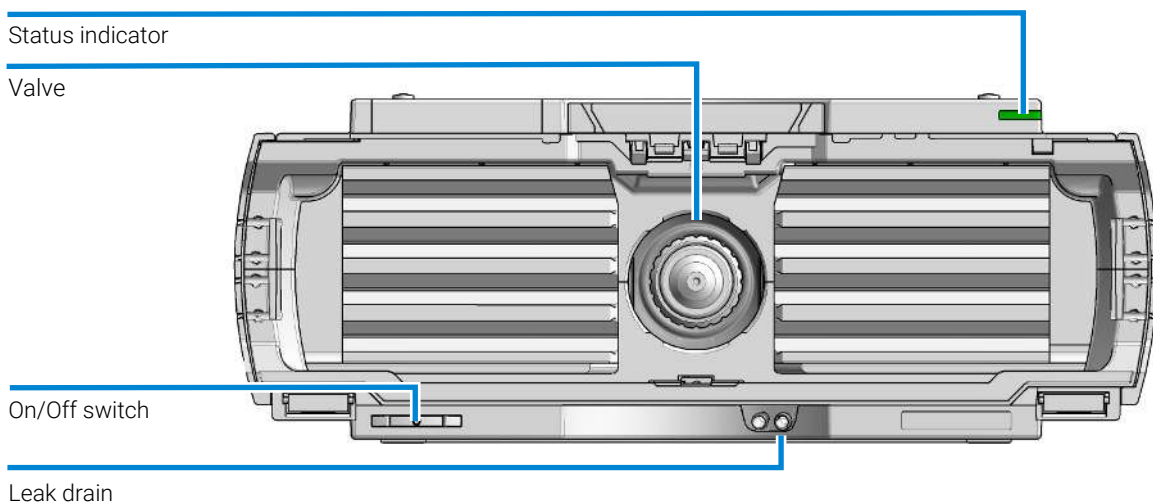
**Figure 5:** Overview of the Vialsampler

## Product Description of the 1290 Infinity III Multicolumn Thermostat (G7116B)

The Agilent 1290 Infinity III Multicolumn Thermostat (MCT) facilitates precise column thermostating over a broad temperature range with cooling down to 20 °C below ambient temperature and heating up to 110 °C.

This capability provides high flexibility for optimized speed and enhanced selectivity of LC separations. Exchangeable ultra-high-pressure valves enable a wide range of applications such as column selection from eight columns in a single MCT, sample preparation for analyte enrichment or matrix removal, alternating column regeneration – and many more.

The MCT matches perfectly with all InfinityLab LC Series systems and can also be combined with previous 1260 and 1290 Infinity Series modules.



**Figure 6:** Overview of the Multicolumn Thermostat

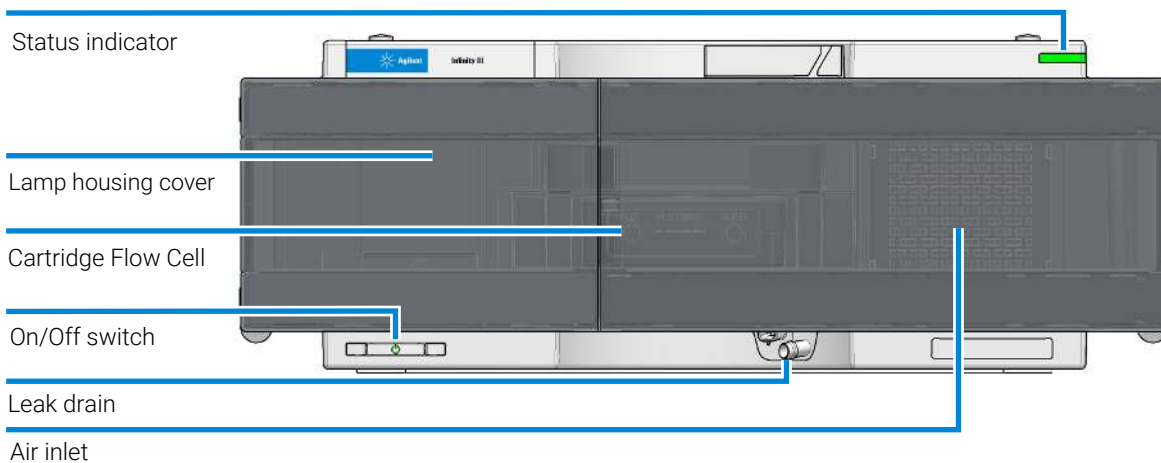
## Product Description of the 1290 Infinity III Diode Array Detector (G7117B)

The Agilent 1290 Infinity III Diode Array Detector (with variable slit) is based on the Agilent Max-Light cartridge cell with optofluidic waveguides that improve light transmission to near 100% efficiency without sacrificing resolution caused by cell dispersions effects.

With typical detector noise levels of  $< \pm 0.6 \mu\text{AU}/\text{cm}$  the 60 mm flow cell gives up to 10 times higher sensitivity than detectors with conventional flow cells.

Any compromising refractive index and thermal effects are almost completely eliminated, resulting in significantly less baseline drift for more reliable and precise peak integration.

For fast separations, this detector has multiple wavelength and full spectral detection at sampling rates up to 240 Hz.



**Figure 7:** Overview of the G7117B Detector

## Product Description of the 1290 Infinity III Variable Wavelength Detector (G7114B)

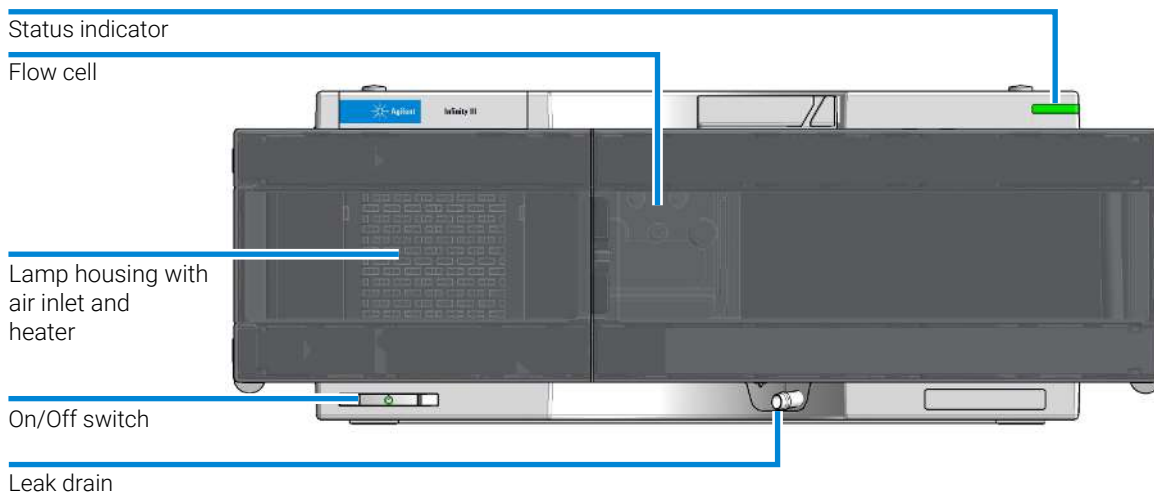
The Agilent 1290 Infinity III Variable Wavelength Detector (VWD) is the most sensitive and fastest detector in its class.

Time-programmable wavelength switching provides sensitivity and selectivity for your applications.

More sample information can be acquired in the dual wavelength mode.

Low detector noise ( $\leq \pm 1.5 \mu\text{AU}$ ) and baseline drift ( $< 1 \cdot 10^{-4} \text{ AU/h}$ ) facilitates precise quantification of trace levels components.

High productivity can be achieved with fast analysis at up to 240 Hz data rates.



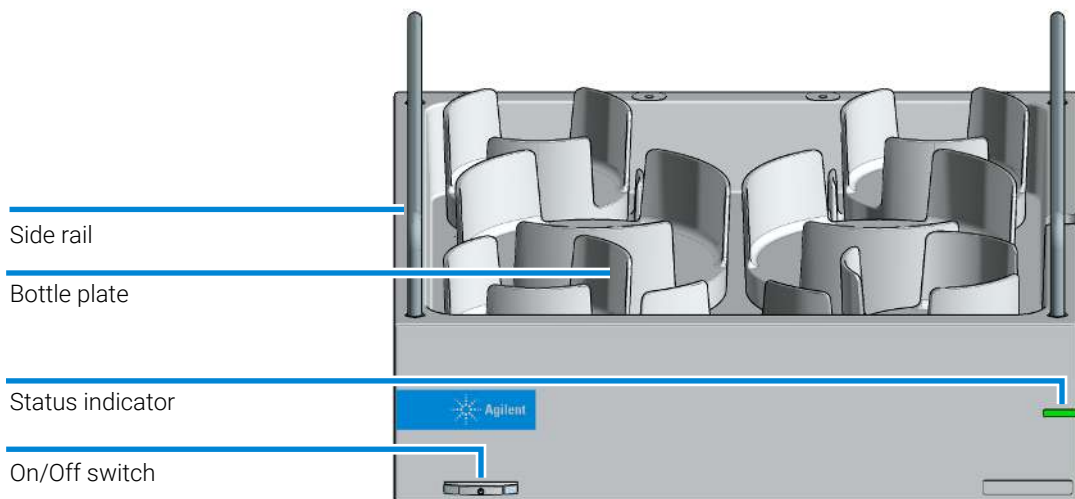
**Figure 8:** Overview of the G7114B Detector



## Product Description of the InfinityLab Level Sensing (G7175A)

The Agilent InfinityLab Level Sensing prevents HPLC instrument or column damage, and time-consuming re-analysis because the mobile phase has run dry. This easy, straightforward system, when used together with OpenLab CDS, also offers a solvent prediction. This feature predicts your HPLC solvent consumption for a complete sequence and warns you before you start so that the injections complete successfully.

The InfinityLab Level Sensing module simplifies the process of LC system preparation and control during analysis. Accurate monitoring of solvent levels ensures optimal productivity.



**Figure 9:** Overview of the Level Sensing module

## Solutions

### 2D-LC Solution

The Agilent InfinityLab 2D-LC Solutions matches separation performance with sample complexity. The solution offers intuitive software, completely pre-configured systems, and an easy starter kit.

Furthermore an easy switch between 1D-UHPLC and 2D-LC guarantees you highest usage of a single system.

The software is designed for fastest method setup in all available modes, whether it is heart-cutting 2D-LC, multiple heart-cutting 2D-LC or comprehensive 2D-LC.

Application areas are numerous and include pharmaceutical and biopharmaceutical (impurity) profiling, all -omics areas, analysis of food matrices, herbal medicine, polymer analysis, flavor analysis and many more.

### High Dynamic Range DAD Solution

The Agilent 1290 Infinity III High Dynamic Range DAD Solution combines the signals from two diode array detectors (DAD). The DADs are assembled with different path length Max-Light flow cells. This combination offers a 30x wider linear UV-range.

The increased linear UV-range allows detection and quantification of all sample components in a single run, thus making it ideal for analysis of mixtures with widely different concentration levels.

## Method Transfer Solution

The 1290 Infinity III Method Transfer Solution allows you to execute any legacy HPLC or latest UHPLC methods while achieving the same chromatographic results. It provides a seamless method transfer between LCs, regardless of brand.

The Intelligent System Emulation Technology (ISET) emulates the LC system on which the original method was developed, thus achieving the same retention times and peak resolution. Speed up your method development with UHPLC performance and then fine-tune your method by emulating the target system. Run your legacy methods with ISET while taking full advantage of the UHPLC speed, resolution, and sensitivity of the 1290 Infinity III LC.

## 2

# Installation

This chapter provides information on unpacking, checking on completeness, stack considerations and installation of the module.

### **Installing the System Modules 21**

#### **Optimizing the Stack Configuration 22**

Agilent InfinityLab Flex Bench 23

One Stack Configuration 25

Vialsampler One Stack Configuration 26

Two Stack Configuration 27

Vialsampler Two Stack Configuration 29

### **Integration Into the Network 30**

#### **Standard Capillary Setup 31**

#### **Handling Leak and Waste 33**

Drain Connectors Installation 37

Waste Concept 41

Waste Guidance 41

Leak Sensor 42

Handling Leak and Waste in a Mixed Configuration 43

## Installing the System Modules

For details of installation procedures for the modules, refer to the individual module manuals. These manuals also contain information on specifications, maintenance and parts.

## Optimizing the Stack Configuration

You can ensure optimum performance by installing the system in one of the following configurations. These configurations optimize the system flow path, ensuring minimum delay volume.

The following configurations are possible:

- InfinityLab Flex Bench
- Single Stack (maximal 4 modules, in a bench rack or directly on the bench)
- InfinityLab Benchtop, providing more flexibility
- Two Stacks (in a bench rack or directly on the bench)

The table below summarizes the advantages of the different prescribed configurations.

**Table 1:** Overview on pros and cons of different stack configurations

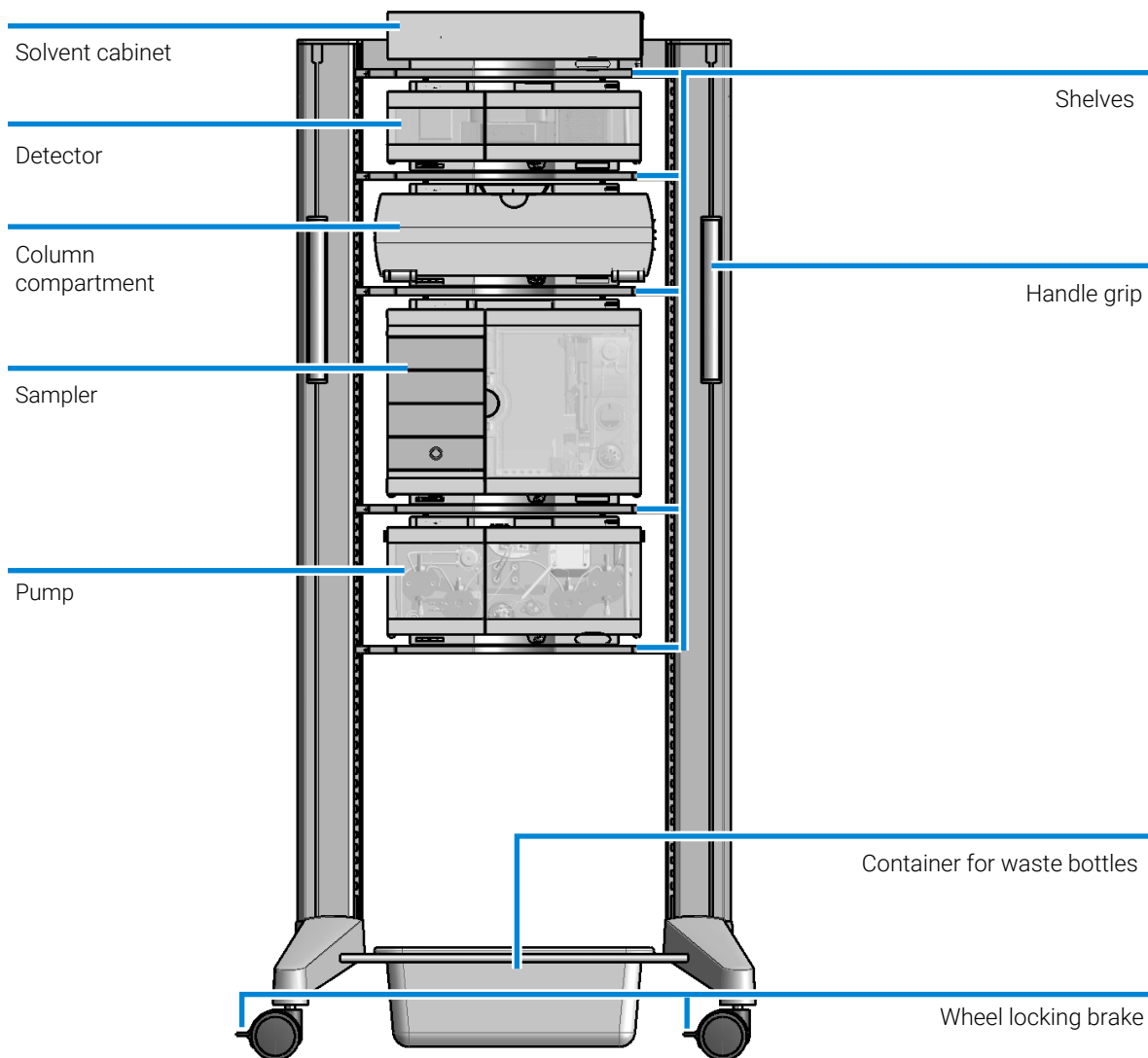
Modules in a stack	InfinityLab Flex Bench Configuration	Single Stack Configuration	Two Stacks Configuration
fewer than 5	Pros <ul style="list-style-type: none"><li>• no bench required</li><li>• mobile</li><li>• optimal access to the modules, solvent bottles, pumps, columns, and accessories</li><li>• integrated waste concept</li></ul>	Pros <ul style="list-style-type: none"><li>• minimal bench space required</li></ul> Cons <ul style="list-style-type: none"><li>• high stack</li></ul>	Pros <ul style="list-style-type: none"><li>• lower stacks</li><li>• flexible combinations</li></ul> Cons <ul style="list-style-type: none"><li>• maximum bench space required</li></ul>
5 and more	+ possible	- not possible	+ possible

## **Agilent InfinityLab Flex Bench**

Agilent recommends using the InfinityLab Flex Bench for all Agilent LC systems.

Main features:

- Increases flexibility in the lab
- Safe moving of LC
- Easy stack customization
- Included waste management

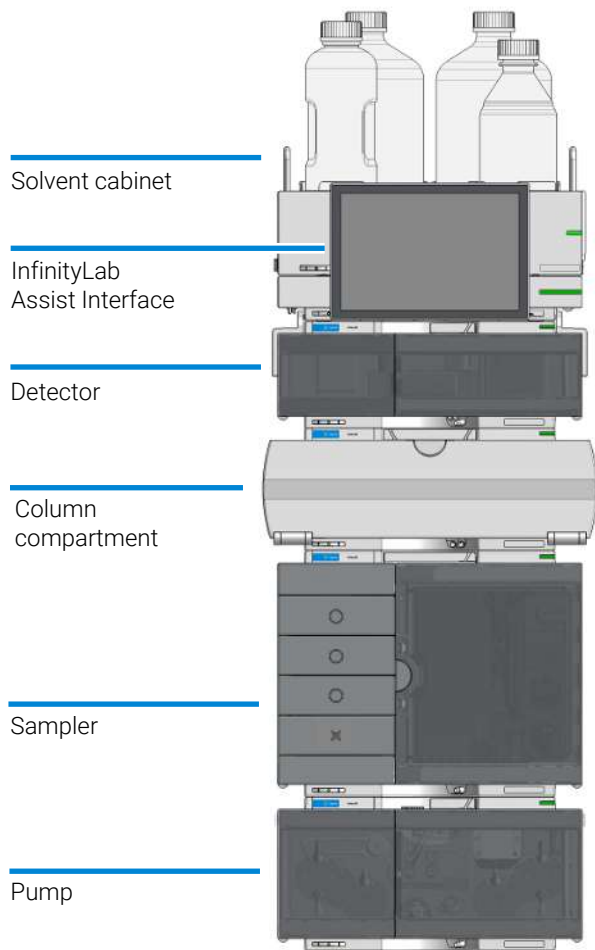


**Figure 10:** Agilent InfinityLab Flex Bench



## One Stack Configuration

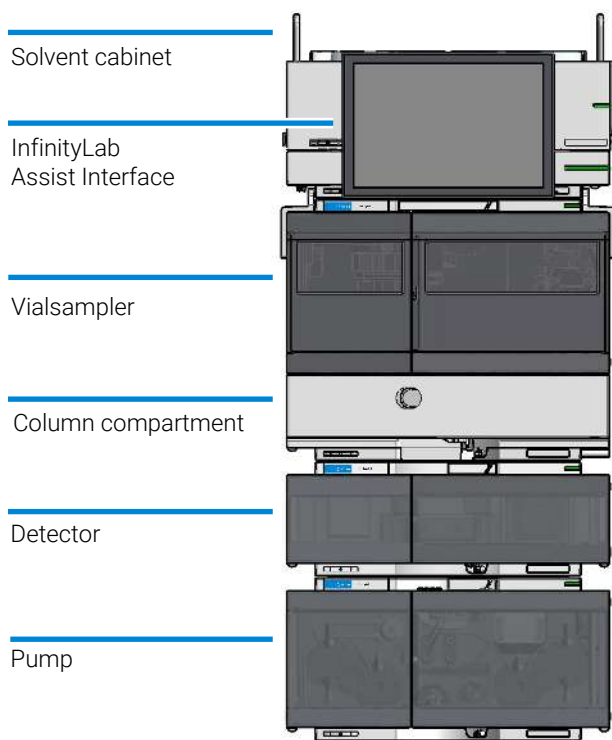
Ensure optimum performance by stacking the modules as shown exemplarily in [Figure 11](#) on page 25. This configuration optimizes the flow path for minimum delay volume and minimizes the bench space required.



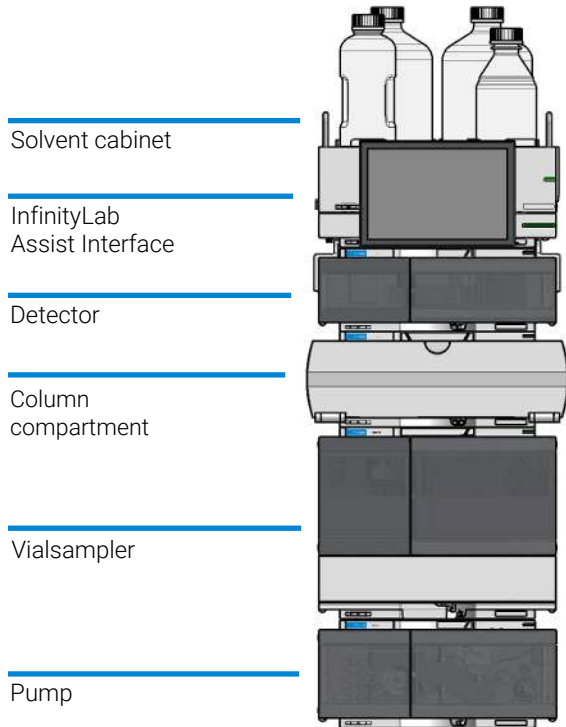
**Figure 11:** Single stack configuration (bench installation, example shows a Multisampler)

## Vialsampler One Stack Configuration

Ensure optimum performance by stacking the modules as shown exemplarily in [Figure 12](#) on page 26 and [Figure 13](#) on page 27 for a Vialsampler set up with and without an Integrated Column Compartment (ICC), respectively. This configuration optimizes the flow path for minimum delay volume and minimizes the bench space required.



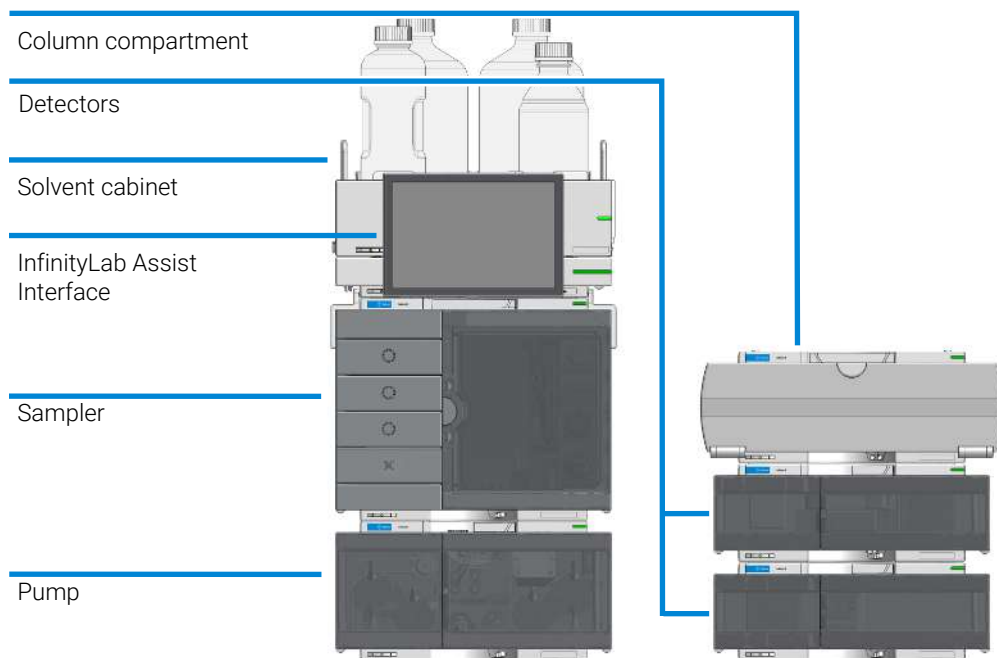
**Figure 12:** Single stack configuration (bench installation, example shows a vialsampler with optional ICC installed)



**Figure 13:** Single stack configuration (bench installation, example shows a Viialsampler coupled with a standalone column compartment)

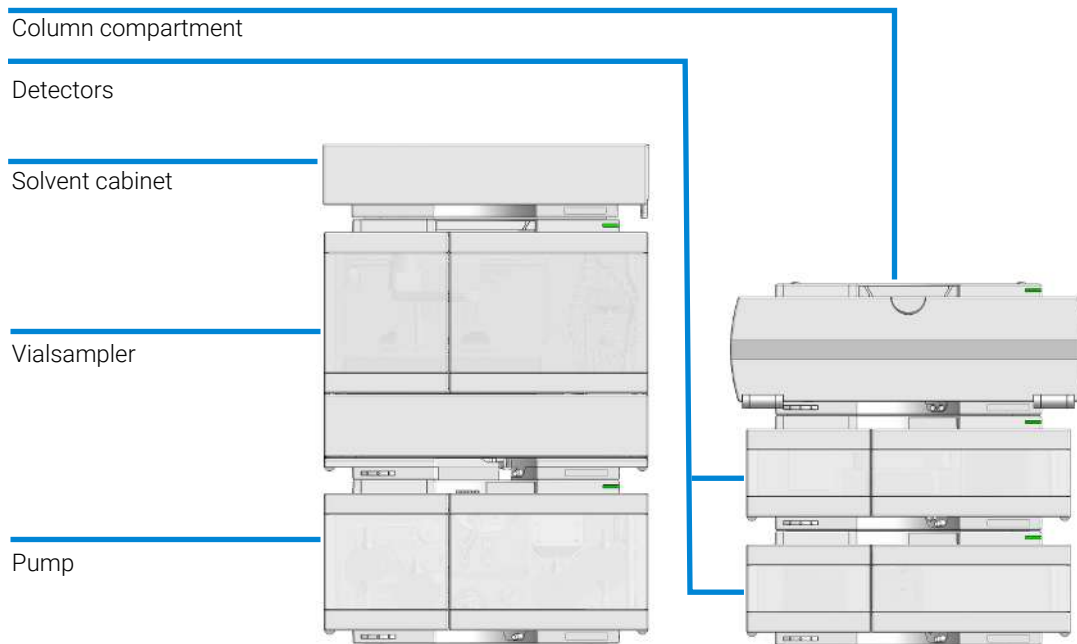
## Two Stack Configuration

To avoid excessive height of the stack (for example when using the system in combination with an additional detector), it is recommended to form two stacks.



**Figure 14:** Two stack configuration (bench installation, example shows a Multisampler)

## Vialsampler Two Stack Configuration

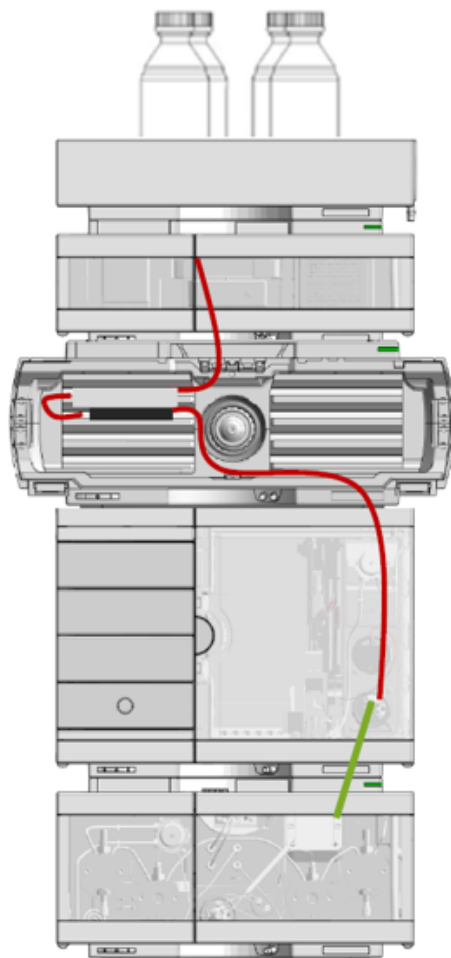


**Figure 15:** Two stack configuration (bench installation, example shows a Vialsampler coupled with a standalone column compartment)

## Integration Into the Network

For network integration of your system refer to user manuals of your modules (chapter *LAN Configuration* ).

## Standard Capillary Setup



**Figure 16:** Standard capillary setup

## Installation

### Standard Capillary Setup

**Table 2:** Capillary properties

Color	Inner diameter	μL per cm
Red	0.12 mm (0.005 inch)	0.113
Green	0.17 mm (0.007 inch)	0.227
Blue	0.25 mm (0.01 inch)	0.491
Orange	0.50 mm (0.02 inch)	



## Handling Leak and Waste

The Agilent InfinityLab LC Series has been designed for safe leak and waste handling. It is important that all security concepts are understood and instructions are carefully followed.

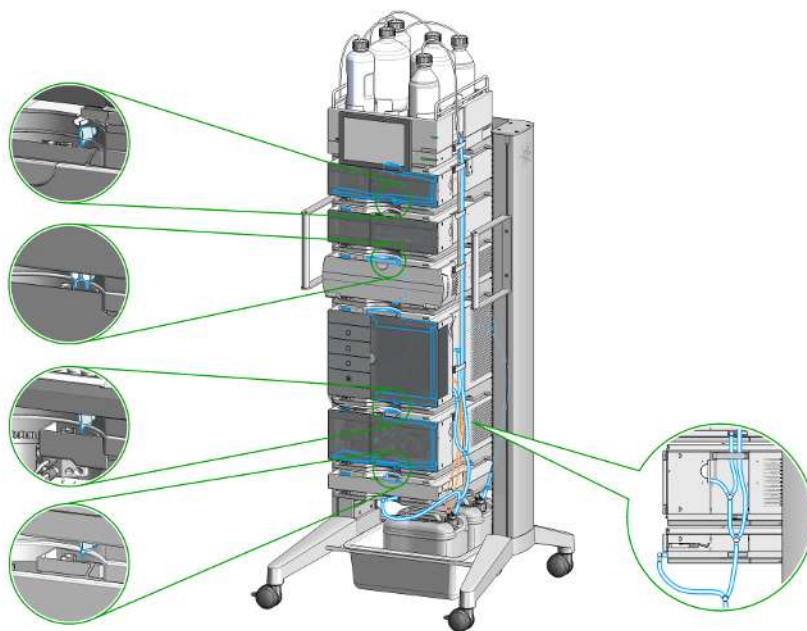
The solvent cabinet is designed to store a maximum volume of 8 L solvent. The maximum volume for an individual bottle stored in the solvent cabinet should not exceed 2 L. For details, see the usage guideline for the Agilent Infinity III Solvent Cabinets (a printed copy of the guideline has been shipped with the solvent cabinet, electronic copies are available on the Internet).

All leak plane outlets are situated in a consistent position so that all Infinity and Infinity II/III modules can be stacked on top of each other. Waste tubes are guided through a channel on the right hand side of the instrument, keeping the front access clear from tubes.

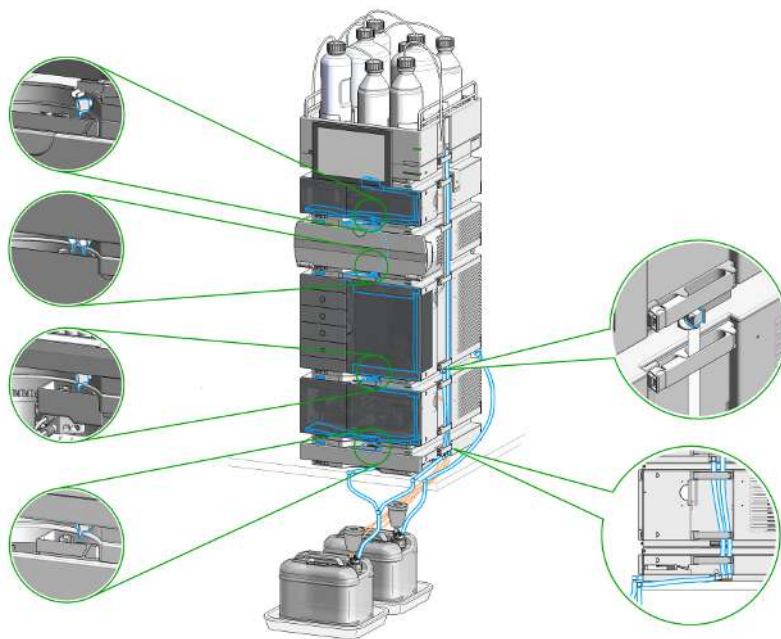
The leak plane provides leak management by catching all internal liquid leaks, guiding them to the leak sensor for leak detection, and passing them on to the next module below, if the leak sensor fails. The leak sensor in the leak plane stops the running system as soon as the leak detection level is reached.

Solvent and condensate is guided through the waste channel into the waste container:

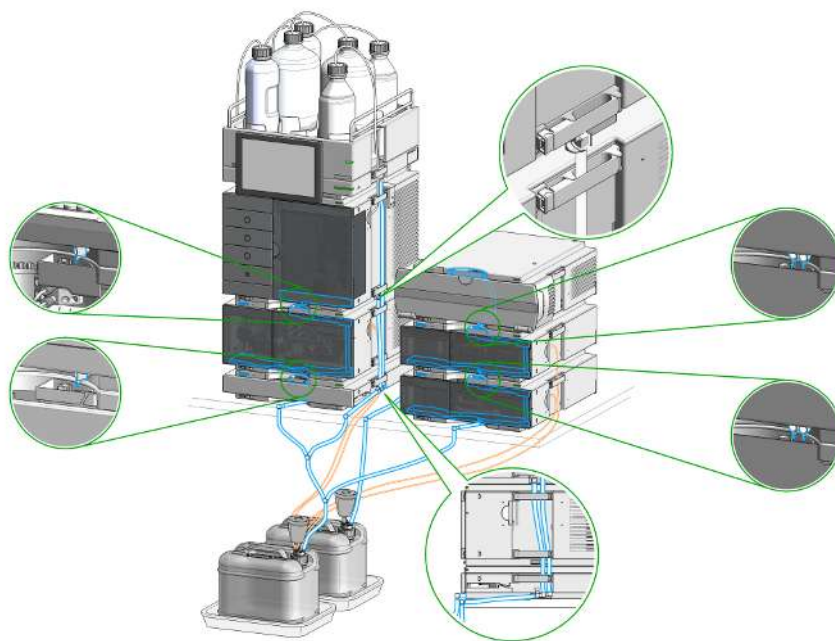
- from the detector's flow cell outlet
- from the Multisampler needle wash port
- from the Sample Thermostat (condensate)
- from the pump's Seal Wash Sensor (if applicable)
- from the pump's Purge Valve or Multipurpose Valve



**Figure 17:** Infinity III Leak Waste Concept (Flex Bench installation)



**Figure 18:** Infinity III Single Stack Leak Waste Concept (bench installation)



**Figure 19:** Infinity III Two Stack Leak Waste Concept (bench installation)

The waste tube connected to the leak plane outlet on each of the bottom instruments guides the solvent to a suitable waste container.

Drain Connectors Installation

Drain Connectors have been developed to improve leak drainage for low flow leaks of high viscosity solvents (for example, isopropanol) in Agilent InfinityLab LC Series Systems. Install these parts to modules where they are missing (usually preinstalled).

- Make sure that dripping adapters are correctly installed on each module in the LC stack, excluding lowest module.
- Remove the dripping adapter if it is appeared to be installed on the lowest module in the LC stack and connect waste tube instead.
- Consider 5004-0000 (Drain Connectors Kit) if drain adaptor is missing on some module(s).

For illustration, see [Handling Leak and Waste](#) on page 33.

Parts required

Qty.	p/n	Description
1	 5004-0000	Drain Connectors Kit

Content of Drain Connectors Kit (p/n 5004-0000)

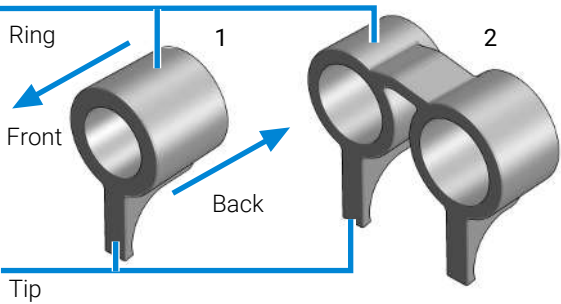




Figure 20: Overview of Drain Connectors: Single (left) and Double (right)

#	Qty.	p/n	Description
Parts can be ordered only as a complete kit.			
1	3	 5043-1834	Single Drain Connector ID3.0-Long
2	1	 5043-1836	Double Drain Connector-Long

**Table 3:** Compatibility of drain connectors and modules

Drain Connector Type	Compatible Module	Compatible Module Type
Double	G7116A/B	Column Compartment
Single	G7114A/B	Detector
	G7115A	
	G7117A/B/C	
	G7121A/B	
	G7162A/B	
	G7165A	
	G7129A/B/C	Sampler
	G7167A/B/C	
	G5668A	
	G7137A	
	G7157A	
	G4767A	
	G7122A	Degasser
	G7104A/C	
	G7110B	Pump
	G7111A/B	
	G7112B	
	G7120A	
	G7131A/C	
	G7132A	
	G5654A	
	G4782A	

**Prerequisites**

- Leak drains of LC modules are clean and free of salt or solvent residuals.

**NOTE**

Do not install drain connectors on the bottom modules of the stack. Drain outlet of the bottom module has to be connected via waste tubing to a suitable waste container (see Leak and Waste Handling in the manual for a respective module).

**NOTE**

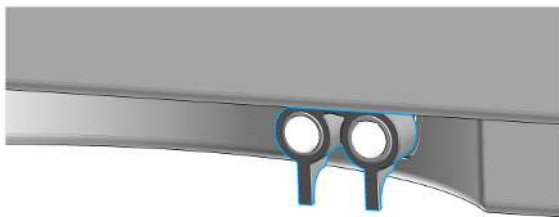
In case of incorrect installation, drain connectors cannot fully perform the intended function.

**NOTE**

It is not required to power off the HPLC stack to install Single and Double Drain Connectors. The installation of the connectors does not affect the analysis performed during the installation.

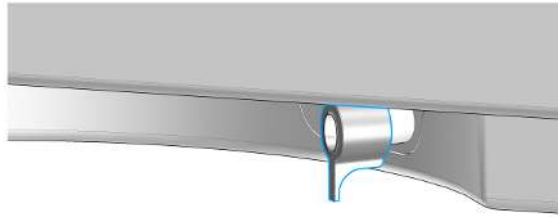
**Install the Double Drain Connector on the leak drain of the  
1260 Infinity III Multicolumn Thermostat (G7116A)/  
1290 Infinity III Multicolumn Thermostat (G7116B)**

- 1 Align the rings with the leak drain outlets of the module, press slightly with the fingers, and slide the connector along the leak drain outlets until it is aligned with the front of the leak drain.



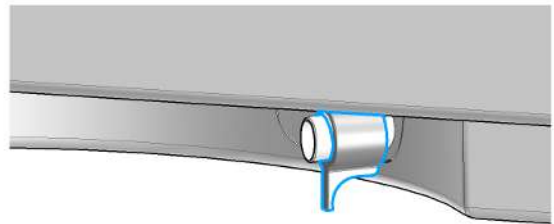
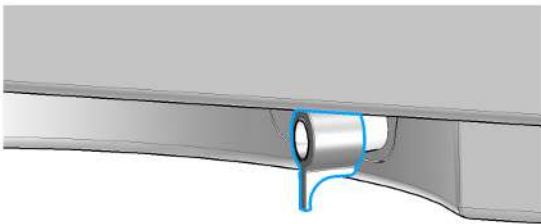
**Install Single Drain Connectors on other modules in the LC stack**

- 1 Align the ring with the leak drain outlet of the module, press slightly with the fingers, and slide the connector along the leak drain outlet until it is aligned with the front of the leak drain.



Make sure that the following requirements are covered:

- The tip of the drain connector points straight down.
- The leak drain outlets and the drain connectors are aligned properly.





## Waste Concept

Agilent recommends using the 5043-1221 (6 L waste can with 1 Stay Safe cap GL45 with 4 ports) for optimal and safe waste disposal. If you decide to use your own waste solution, make sure that the tubes don't immerse in the liquid.



## Waste Guidance



**NOTE**

The waste drainage must go straight into the waste containers. The waste flow must not be restricted at bends or joints.

## Leak Sensor

**CAUTION****Solvent incompatibility**

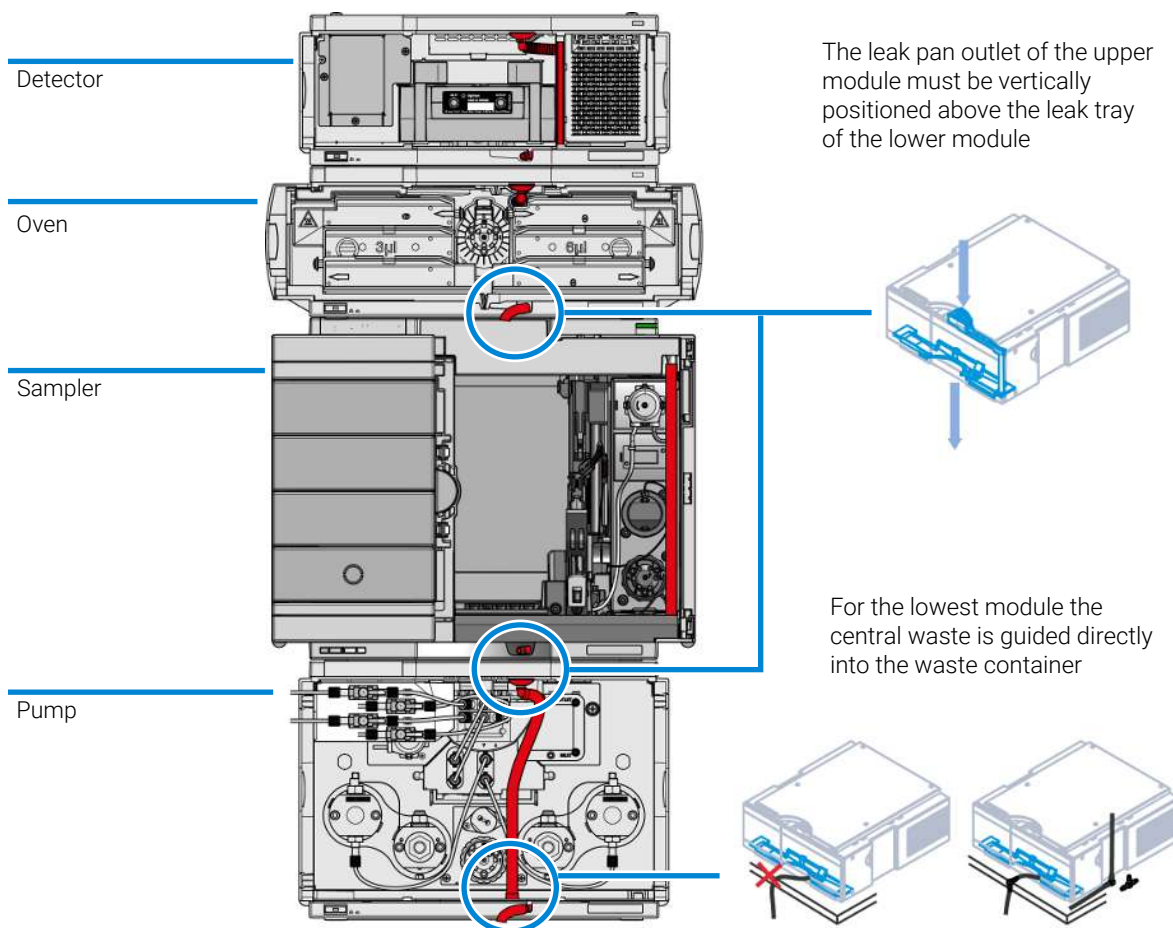
The solvent DMF (dimethylformamide) leads to corrosion of the leak sensor. The material of the leak sensor, PVDF (polyvinylidene fluoride), is incompatible with DMF.

- Do not use DMF as mobile phase.
- Check the leak sensor regularly for corrosion.

**NOTE**

The leak sensor in the sampler is hidden under the ICC Column Heater or Column Shelf respectively.

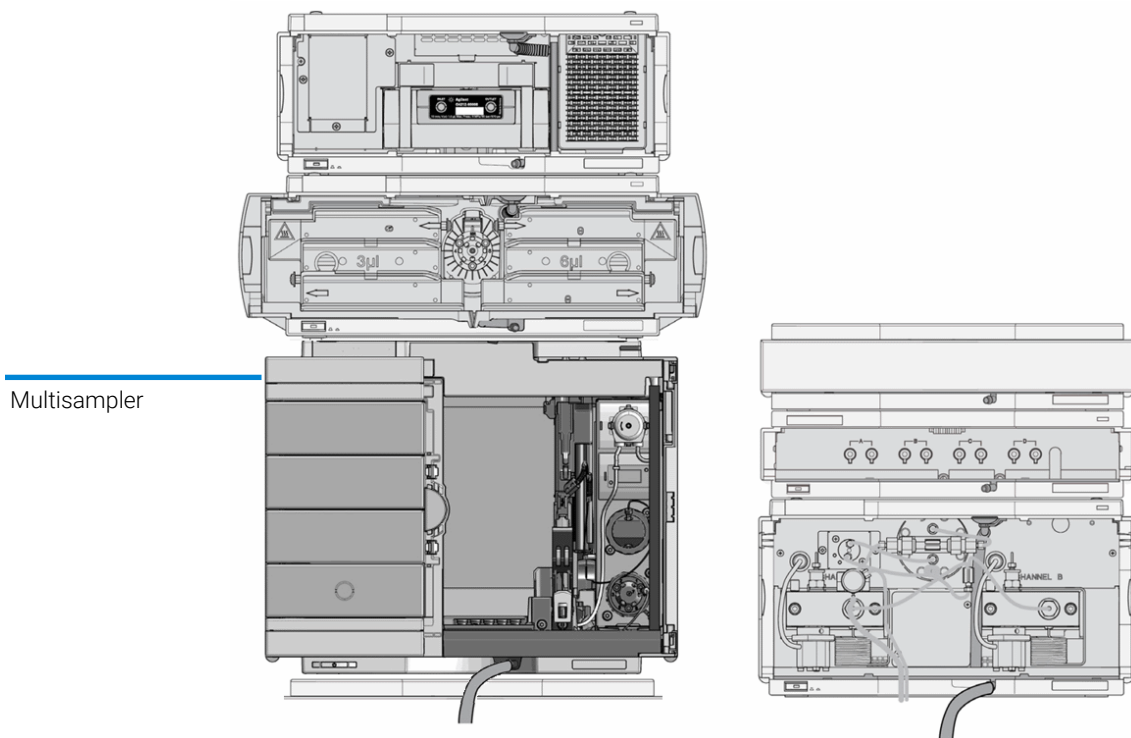
## Handling Leak and Waste in a Mixed Configuration



**Figure 21:** Leak and waste handling with multisampler in a mixed configuration as an example

### NOTE

Flush solvent from the washport of the multisampler is guided out to the right of the instrument.



**Figure 22:** Leak and waste handling with multisampler in a mixed configuration as an example (two stack configuration)

#### NOTE

Do not place the multisampler directly on the bench if a sample cooler or sample thermostat is installed.



## 3 Configuration Settings

This chapter describes how to configure the system.

**General Information on LAN Configuration 46**

**Instrument Configuration 47**

**Lab Advisor 49**

Adding a New System 49

Installing Add-ons 52

## General Information on LAN Configuration

LAN configuration is executed from the module with direct LAN connection to the controller software. This must be the module (usually the detector) with the highest data rate.

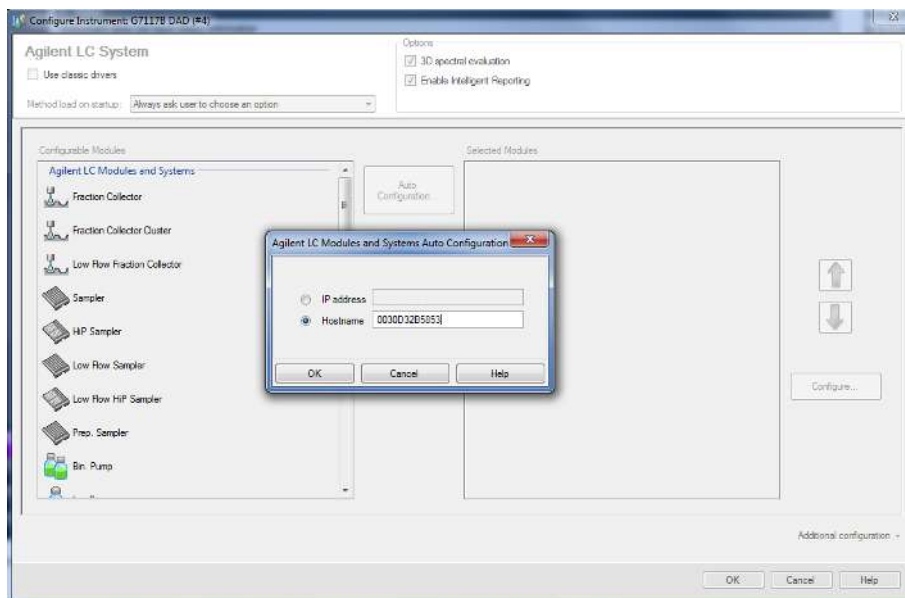
## Instrument Configuration

Example shows an instrument configuration with a Diode Array Detector.

- 1 Set the switches of the Configuration switch at the rear of the module:
  - a All switches DOWN: module uses the default IP address 192.168.254.11.



- b Switch 4 UP and others DOWN: module uses DHCP.
  - c Switch 5 UP and others DOWN: modules uses STORED address.
- 2 Enter the setup information (MAC <sup>1</sup> / IP address and/or Instrument Name).
  - a Agilent OpenLab ChemStation (Configure Instrument):



<sup>1</sup> MAC address can only be used in DHCP DIP-switch configuration.

## Configuration Settings

### Instrument Configuration


b Lab Advisor (Instrument Overview - Add Instrument):

**System Properties**

**System**

System Name:  Description:

**Instruments**

Instrument Name:	Instrument Address:	Instrument Type:
 <input type="text" value="G7117B"/>	<input type="text" value="0030D32B5853"/> <input type="text" value="192.168.254.11"/>	<input type="text" value="Agilent LC/CE"/>

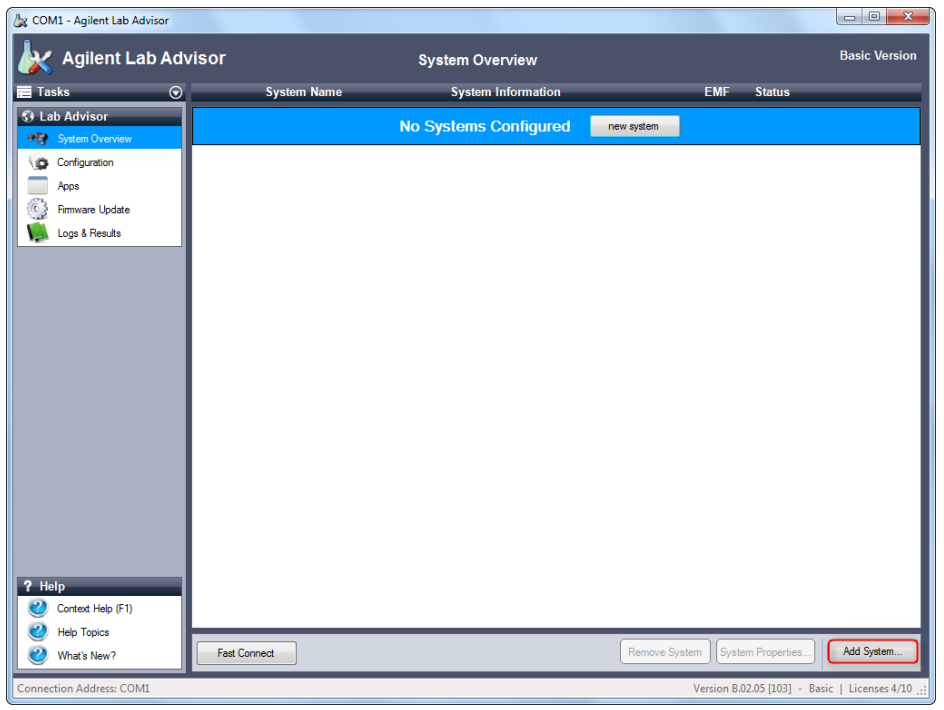
Reconnect: ☐



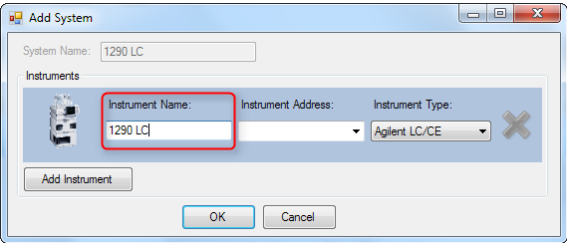
# Lab Advisor

## Adding a New System

- 1 In the Action Panel of the System Overview, click Add System.



The Add System dialog box is displayed.

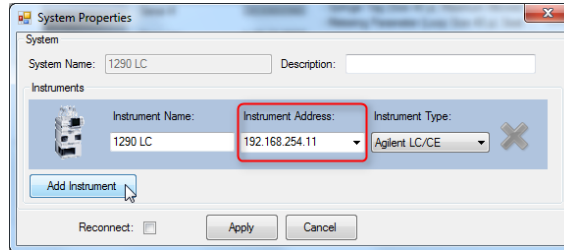


- 2 Enter a name in the **Instrument Name** field.

#### NOTE

If your system comprises just one instrument, the **Instrument Name** is copied to the **System Name** field.

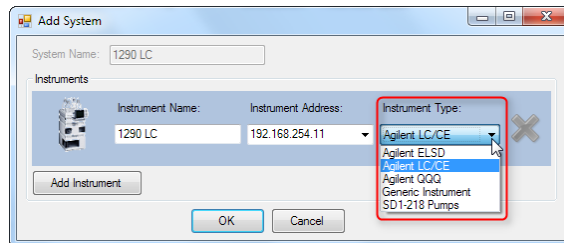
- 3 Enter the connection details in the **Instrument Address** field.



#### NOTE

The **Instrument Address** can be an IP address, the host name or, if you are connecting using a serial cable, the COM port.

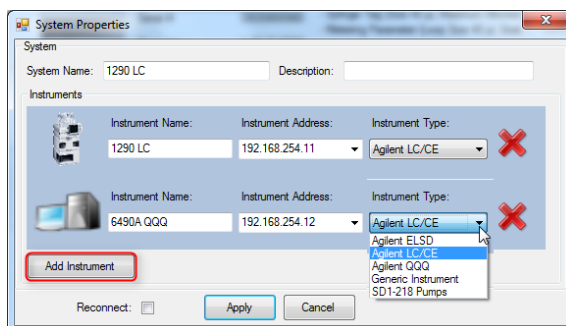
- 4 Click the **Instrument Type** down-arrow and select the type of instrument you are adding from the list. The default setting is **Agilent LC/CE**. Additional instrument types become available when the respective add-ons are installed.



#### NOTE

By default, the **Instrument Type** drop-down list contains only the entry **Agilent LC/CE**. Additional instrument types can be added by installing the respective add-ons (see [Installing Add-ons](#) on page 52).

- 5 If your system comprises more than one instrument, click **Add Instrument** and complete the details as above.

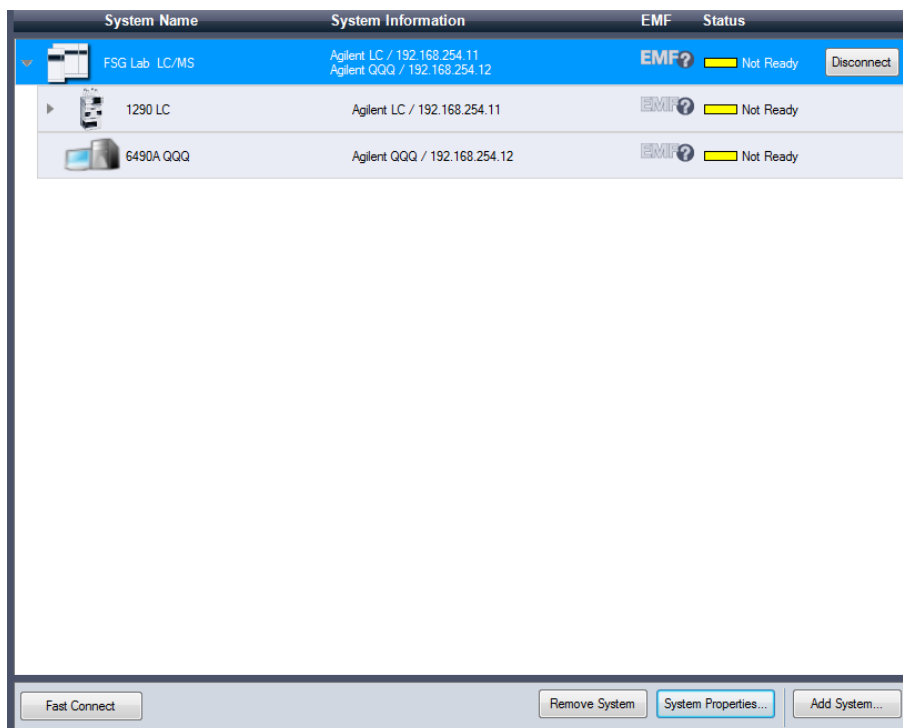


## NOTE

As soon as you add a second instrument, the **System Name** field is activated to allow you to edit the system name.

- 6 Click **OK** to finish adding the system and close the **Add System** dialog box.

The system becomes visible in the **System Overview**, and Lab Advisor tries to connect to it.



## Installing Add-ons

Add-ons are installed from the **Configuration** screen, using a Lab Advisor Extension file with the extension .LAX.

### NOTE

You need Administrator rights in order to install Add-ons.

- 1 In the Global Tasks section of the Navigation Panel, click **Configuration**.  
The **Configuration** screen is displayed.
- 2 Click **Add-ons** to navigate to the **Configuration - Add-ons** screen.

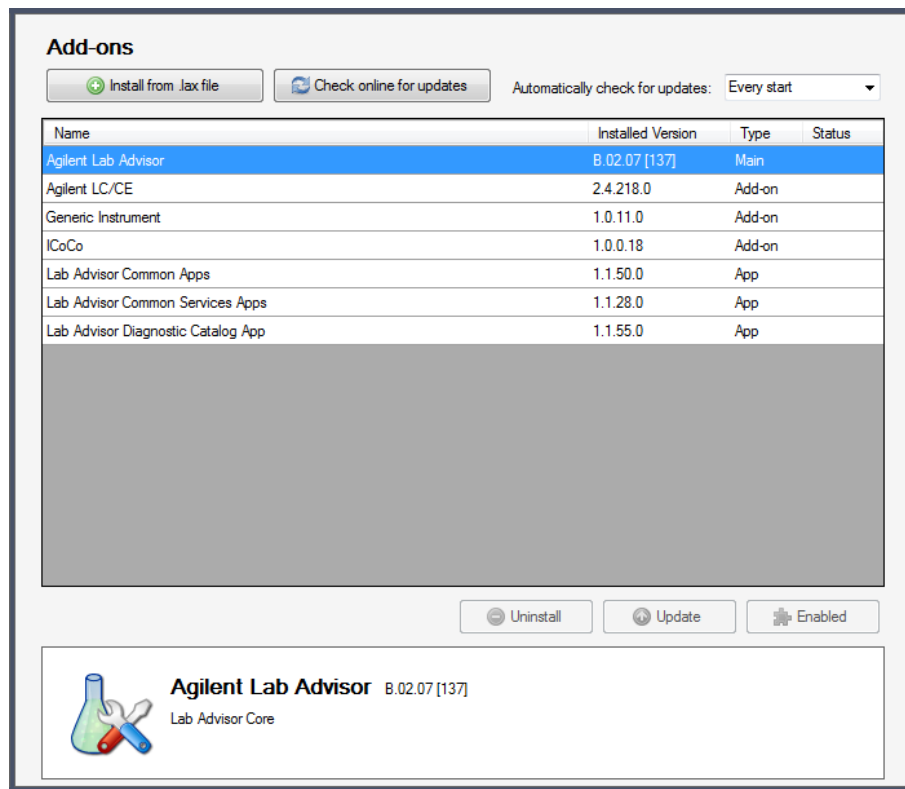


Figure 23: Add-ons in Configuration

The **Configuration - Add-ons** screen contains a table listing all the Add-ons that are already installed.

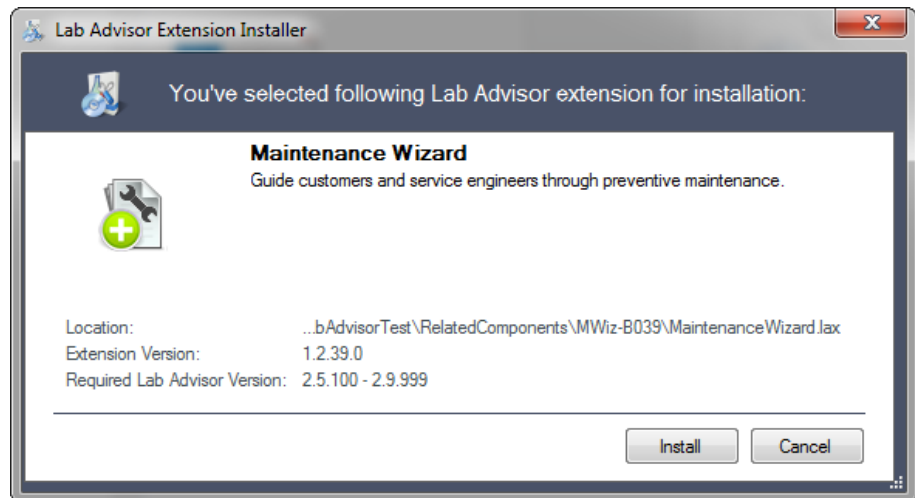
**3** Click **Install from .lax file**.

A file selection dialog box is displayed to allow you to select the App or Add-on to install.

**4** Navigate to the folder containing the Add-on files, select the .lax file and click **Open** to install the Add-on.

**5** Click **Yes** when the request to shut down Lab Advisor appears.

Lab Advisor shuts down and the Add-on installation is started.



When the installation is finished, the newly installed Add-on is included in the table in the **Configuration - Add-ons** screen.

# 4

## Quick Start Guide

This chapter provides information on running an Agilent 1290 Infinity III LC System.

### **Best Practices 55**

Daily / Weekly Tasks 55

Power Up / Shut Down the Pump 55

Prepare the Pump 56

How to Deal With Solvents 57

Select Channels for Gradient Valve (MCGV/4CGV) 57

Optional Inline Filter 57

### **Prepare a Run 59**

### **Installation Checkout (Customer presence recommended) 69**

Checkout Columns 70

Checkout Samples 71

Checkout Method for Isocratic, 400 bar and FLD only systems 73

Checkout Method for 1260 and 1290 Systems with UV Detectors 76

Checkout Method for 1260 and 1290 Systems with ELSD or RID 80

## Best Practices

### Daily / Weekly Tasks

#### Daily Tasks

- Replace solvents and solvent bottles for mobile phases based on water/ buffer.
- Replace solvents and solvent bottles for organic mobile phase latest every second day.
- Check presence of seal wash solvent.
- Purge each channel with fresh solvent at 2.5 – 3 mL/min for 5 min before operation.
- Equilibrate the system with composition and flow rate of subsequent method.

#### Weekly Tasks

- Change seal wash solvent (10 % isopropanol in water) and bottle.
- If applications with salts were used, flush all channels with water and remove possible salt deposits manually.
- Inspect solvent filters for dirt or blockages. Exchange if no flow is coming out of the solvent line when removed from the degasser inlet.

### Power Up / Shut Down the Pump

#### Prepare the Pump

- Use fresh or different mobile phase (as required).
- Purge each channel with 2.5 – 3 mL/min for 5 min. Open the manual purge valve or use the purge command, depending on the pump type.

## Long-Term Shut-Down of the System

- Flush system with water to remove buffer.
- Remove all samples from the sampler and store according to good laboratory practice.
- Use recommended solvents to store the system.
- Power off the system.

## Prepare the Pump

### Purge

Use the Purge function to:

- fill the pump,
- exchange a solvent,
- remove air bubbles in tubes and pump heads.

### Condition

Use the Conditioning function:

- daily when starting the pump,
- to minimize pressure ripple by dissolving air bubbles in pump heads.

#### NOTE

Condition your complete system with solvents and composition of your application (for example 50 %/50 % A/B at 0.5 mL/min).

### Seal Wash

Seal Wash guarantees a maximum seal life time. Use Seal Wash:

- When using buffers with elevated salt concentrations
- When using volatile solvents with non-volatile additives



**CAUTION**

Contaminated seal wash solvent

- Do not recycle seal wash solvent to avoid contamination.
- Weekly exchange seal wash solvent.

## How to Deal With Solvents

- Use clean bottles only.
- Exchange water-based solvents daily.
- Select solvent volume to be used up within 1 – 2 days.
- Use only HPLC-grade solvents and water filtered through 0.2 µm filters.
- Label bottles correctly with bottle content, and filling date / expiry date.
- Use solvent inlet filters.
- Reduce risk of algae growth: use brown bottles for aqueous solvents, avoid direct sunlight.

## Select Channels for Gradient Valve (MCGV/4CGV)

- Use lower channels (A and/or D) for buffer solutions.
- Regularly flush all MCGV/4CGV channels with water to remove possible salt deposits.
- Check compatibility of buffers and organic solvents to avoid precipitation.

## Optional Inline Filter

The pump can be equipped with an additional inline filter (5067-5407 (Inline filter assembly, material: stainless steel) or 5720-0003 (Bio Inline Filter for Infinity II Bio Flexible Pumps) ) with a nominal filter pore size of 0.3 µL.

Advantages of the inline filter:

- Very small internal volume

- Specified for working at high pressures
- Possibility of back-flushing the filter

Using the inline filter is recommended:

- to protect the downstream system from blockages,
- for solvent combinations that can form precipitation after mixing,
- for applications running with buffers.

General hints for effective use of the inline filter:

- filter solvents before use,
- follow best practices,
- back-flush the filter weekly,
- exchange the filter frit regularly.

#### CAUTION

##### Damage to the valve

- Use the filter flush mode only if the optional inline filter is installed.

See *1290 Infinity II Inline Filters Technical Note (G7167-Inline-Filters-TechPu-en-SD-29000397.pdf, SD-29000397)* for further reference.

## Prepare a Run

This procedure exemplarily shows how to prepare a run. Parameters as shown in the screenshots may vary, depending on the system installed.

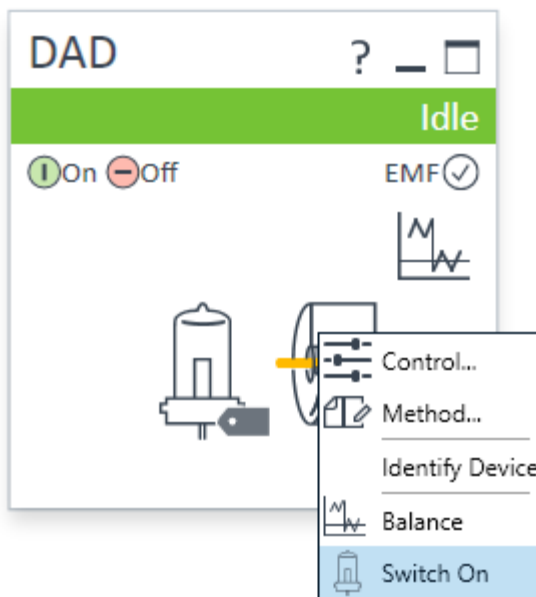
### WARNING

#### Toxic, flammable and hazardous solvents, samples and reagents

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

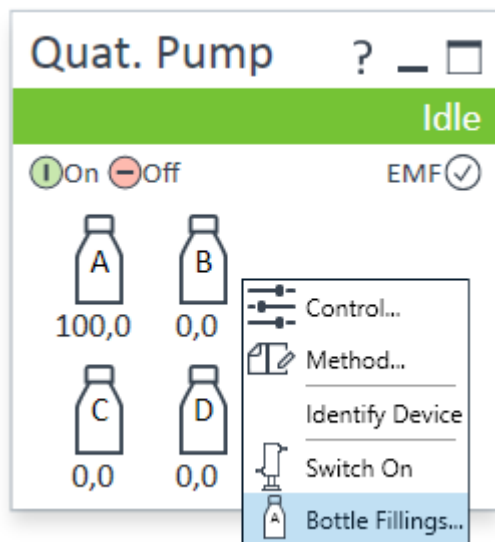
- When working with these substances observe appropriate safety procedures (for example by wearing goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the vendor, and follow good laboratory practice.
- Do not use solvents with an auto-ignition temperature below 200 °C (392 °F). Do not use solvents with a boiling point below 56 °C (133 °F).
- Avoid high vapor concentrations. Keep the solvent temperature at least 40 °C (72 °F) below the boiling point of the solvent used. This includes the solvent temperature in the sample compartment. For the solvents methanol and ethanol keep the solvent temperature at least 25 °C (45 °F) below the boiling point.
- 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.
- Do not use bottles that exceed the maximum permissible volume (2.5 L) as specified in the usage guidelines.
- 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 Switch on the detector.



- 2 Fill the solvent bottles with adequate solvents for your application.
- 3 Place solvent tubings with bottle head assemblies into the solvent bottles.
- 4 Place solvent bottles into the solvent cabinet.

- 5 Solvent bottle filling dialog (in the software).



#### Solvent Bottle

##### Fillings

	Actual Volume		Total Volume	
A:	<input type="text" value="0.80"/>	liter	<input type="text" value="1.00"/>	liter
B:	<input type="text" value="0.92"/>	liter	<input type="text" value="1.00"/>	liter
C:	<input type="text" value="0.78"/>	liter	<input type="text" value="1.00"/>	liter
D:	<input type="text" value="0.81"/>	liter	<input type="text" value="1.00"/>	liter

##### Actions

- ☒ Prevent analysis if level falls below  liter
- ☐ Turn pump off if running out of solvent

Ok

Cancel

Help

- 6 Purge the pump (in normal usage scenario).

**OR:** Prime the pump (after installation of the system).

## NOTE

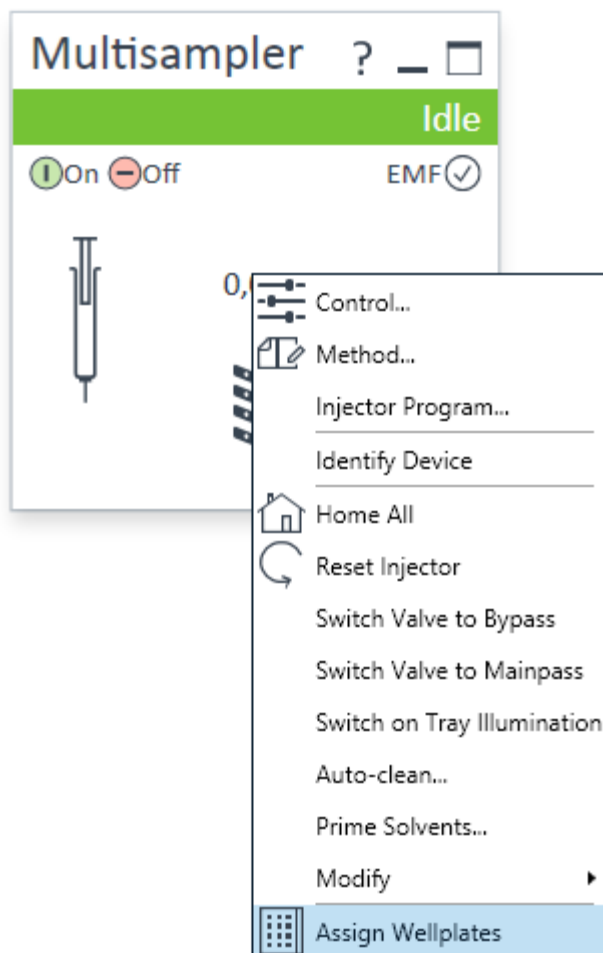
For details on priming and purging, refer to the technical note *Best Practices for Using an Agilent LC System Technical Note (InfinityLab-BestPractice-en-SD-29000194.pdf, SD-29000194)*.

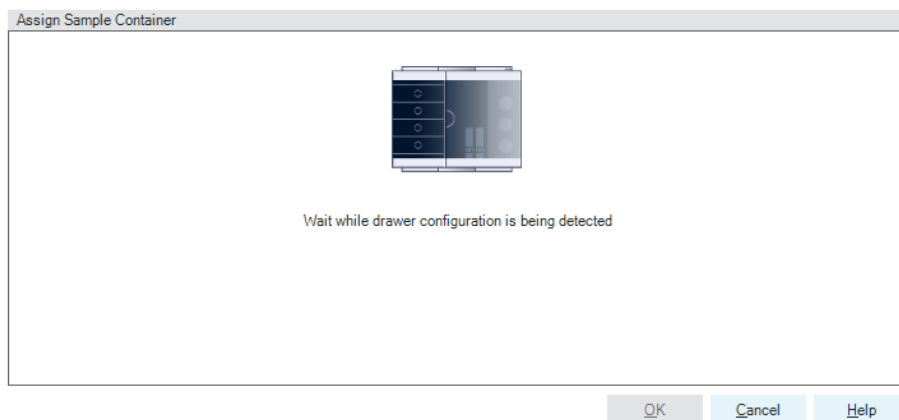
7 Change solvent (if necessary).

The screenshot displays the 'Quat. Pump (G7111B)' software interface. At the top, there are tabs for 'Quat. Pump', 'VWD', 'Sampler', 'DAD', and 'Column Comp.'. The main window is divided into several sections:

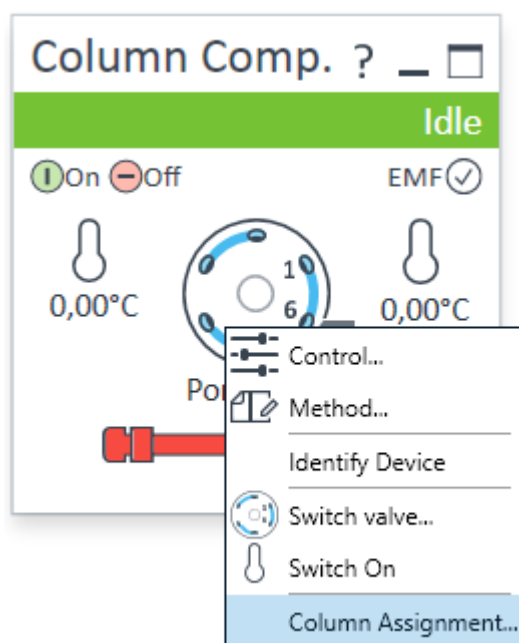
- Flow:** A single input field shows '0.800' mL/min.
- Solvents:** A list of four solvents (A, B, C, D) with their respective percentages and names. Solvent B is selected with a checkmark.
  - A: 60.0 % Water
  - B: 40.0 % Acetonitrile
  - C: 0.0 %
  - D: 0.0 %
- Pressure Limits:** Two input fields for 'Min:' (0.00 bar) and 'Max:' (600.00 bar).
- Stoptime:** A radio button selection between 'As Injector/No Limit' and 'Off'. The 'As Injector/No Limit' option is selected, with a time input of 10.00 min.
- Posttime:** A radio button selection between 'Off' and a time input of 1.00 min. The 'Off' option is selected.
- Advanced:** A sidebar on the right containing:
  - Minimum Stroke:** Radio button selection between 'Automatic' (selected) and a value of 20 µL.
  - Compressibility:** Radio button selection between '95 \*10<sup>-6</sup> / bar' (selected) and 'No compensation'.
  - Maximum Flow Gradient:** An input field showing 100.000 mL/min².
  - Primary Channel:** A dropdown menu set to 'Automatic'.
- Import Timetable...:** A button located below the Solvents section.
- Timetable (1/100 events):** A section at the bottom right with a scrollbar and a play button icon.

- 8 Choose the tray format of the sampler.





- 9 Add a new column.



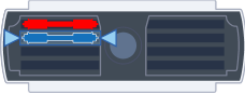


10 Enter the column information.

Plumbing

Valve Position	Location
1	Left 1
2	Left 2

Visualization



Valve Type: 2-pos/6-port valve 600 bar (5067-4137)

Column Tag Information

>>

Location	Color Code	Description	Length [mm]	Diameter [mm]	Particle Size [µm]	Max. Pressure [bar]	Injections
Left 1	Red		0	0,0	0,0	0	0
▶ Left 2	Blue		0	0,0	0,0	0	0
Left 3	None		0	0,0	0,0	0	0
Left 4	None		0	0,0	0,0	0	0
Right 1	None		0	0,0	0,0	0	0
Right 2	None		0	0,0	0,0	0	0
Right 3	None		0	0,0	0,0	0	0
Right 4	None		0	0,0	0,0	0	0

Ok/Write TagCancelHelp

## 11 Select the column position.

Temperature

Left:

☐ Not Controlled  
☒ 40.0 °C  
☐ As Detector Cell  
☐ Unchanged

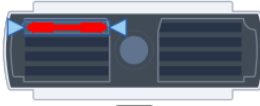

Right:

☐ Not Controlled  
☐ 25.0 °C  
☐ As Detector Cell  
☐ Unchanged  
☒ Combined

Valve Position/Column

☐ Use Current Column / Position  
☒ Use Selected Column / Position

Position 1

☐ Enforce column for run

Stoptime

☒ As Pump/Injector  
☐ 1.00 min

Posttime

☒ Off  
☐ 1.00 min

Advanced

Enable Analysis

☒ when front door open

Left:

☐ With any temperature  
☒ When temperature is within

± 0.8 °C for

0.0 min

Right:

☐ With any temperature  
☒ When temperature is within

± 0.8 °C for

0.0 min

Valve Position/Column After Run

☒ Do not switch  
☐ Switch to position / column at beginning of run  
☐ Increase valve position / column  
☐ Use valve position / column

Position 1

Timetable (empty)

Ok

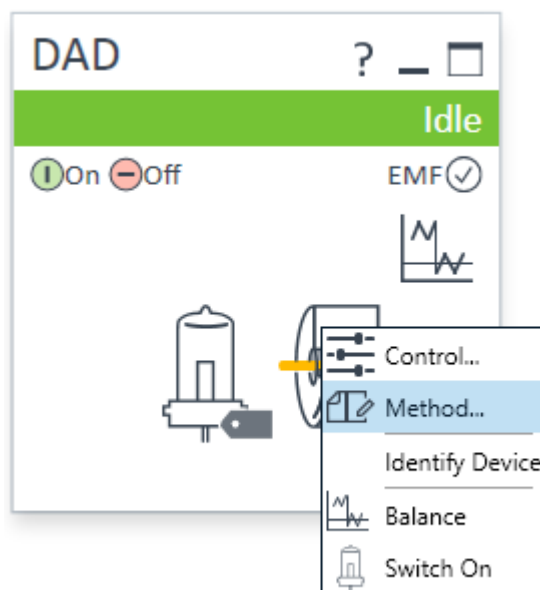
Apply

Cancel


1290 Infinity III LC System Manual

66

12 Set the detector according to the needs of your method.



Quat. Pump | VWD | Sampler | **DAD** | Column Comp.

**DAD (G7115A)** 

**Signals**

	Acquire	Wave length	Band width	Reference Wavelength	Reference Bandwidth
Signal A	<input checked="" type="checkbox"/>	254	4	<input checked="" type="checkbox"/> 360	100 nm
Signal B	<input type="checkbox"/>	254	4	<input type="checkbox"/> 360	100 nm
Signal C	<input type="checkbox"/>	210	4	<input type="checkbox"/> 360	100 nm
Signal D	<input type="checkbox"/>	230	4	<input type="checkbox"/> 360	100 nm
Signal E	<input type="checkbox"/>	280	4	<input type="checkbox"/> 360	100 nm
Signal F	<input type="checkbox"/>	260	4	<input type="checkbox"/> 360	100 nm
Signal G	<input type="checkbox"/>	270	4	<input type="checkbox"/> 360	100 nm
Signal H	<input type="checkbox"/>	290	4	<input type="checkbox"/> 360	100 nm

**Peakwidth**

> 0.0063 min (0.13 s response time) (40 Hz) ▾

**Stoptime** **Posttime**

☒ As Pump/Injector      ☒ Off  
☐ 1.00 min      ☐ 1.00 min

**Advanced**

**Spectrum**

Store: ▾

Range from: 190 to 400 nm

Step: 2.0 nm

**Analog Output**

Zero Offset: 5 %

Attenuation: 1000 mAU

**Margin for negative Absorbance** **Slit**

100 mAU      4 nm

**Autobalance** **Lamps on required for acquisition**

☒ Prerun      ☒ UV Lamp  
☐ Postrun      ☐ Vis Lamp

▶ Timetable (empty)

**NOTE**

For details on running a method, see Set up the Checkout Method as an example.

## Installation Checkout (Customer presence recommended)

For instruments where the checkout columns mentioned in Overview of column options for different pumps are included, the checkout is run after the complete installation of the module stack to affirm the functionality of all modules.

The checkout confirms that each module performs and is connected correctly. The chromatography should show a single peak for FLD, RID or ELSD or four or nine separated peaks, respectively, but is not a substitute for system suitability tests or qualifications.

The checkout for UV and Fluorescent Light Detectors should be run with one of the checkout columns supplied with the pump or with an equivalent column to ensure separation of the compounds.

The checkout for RIDs and ELSDs is done with a 5022-2159 (Restriction capillary, SST 0.12 mm ID, 2 m long) (use 5005-0046 (Capillary MP35N 0.12 mm x 2 m) for 1290 Infinity II Bio LC System).

When Analytical Fraction Collector is installed in the LC System, acetonitrile and water mobile phases with 0.1 % formic acid (as needed in the FC checkout procedure) can be used for the system checkout procedure without any further method modifications.

### NOTE

If a system is equipped with multiple detectors, only run one checkout run and always use the column based procedure for system checkout.

## Checkout Columns

The information in this table applies to Infinity II and Infinity III modules.

**Table 4:** Overview of column options for different pumps

Pump	Option 1	Option 2	Option 3
1260 Isocratic, Quaternary and Binary Pump (G7110B/G7111B/G7112B)	695975-902T, InfinityLab Poroshell 120 EC-C18, 4.6 x 100 mm, 2.7 µm, 600 bar	693975-302T, InfinityLab Poroshell 120 EC-C18, 3.0 x 150 mm, 2.7 µm, 600 bar	699975-302T, InfinityLab Poroshell 120 EC-C18, 3.0 x 50 mm, 2.7 µm, 600 bar
1260 Quaternary Pump VL (G7111A)	695970-902T, InfinityLab Poroshell 120 EC-C18, 4.6 x 100 mm, 4 µm, 600 bar	699975-902T, InfinityLab Poroshell 120 EC-C18, 4.6 x 50 mm, 2.7 µm, 600 bar	693970-902T, InfinityLab Poroshell 120 EC-C18, 4.6 x 150 mm, 4 µm, 600 bar
1260 Flexible Pump (G7104C)	693575-302, InfinityLab Poroshell 120 EC-C18, 3.0 x 150 mm, 2.7 µm, 1000 bar	695575-302, InfinityLab Poroshell 120 EC-C18, 3.0 x 100 mm, 2.7 µm, 1000 bar	699675-902, InfinityLab Poroshell 120 EC-C18, 2.1 x 50 mm, 1.9 µm, 1300 bar
1290 Pumps (G7120A and G7104A)	699675-902, InfinityLab Poroshell 120 EC-C18, 2.1 x 50 mm, 1.9 µm, 1300 bar		
1260 Bio-inert Pump (G5654A)	653750-902, AdvanceBio Peptide Mapping 120Å, 2.1 x 150 mm, 2.7 µm, 600 bar		
1260 and 1290 Bio Flexible Pump and 1290 Bio High-Speed Pump (G7131A/C and G7132A)	691975-302T, InfinityLab Poroshell 120 EC-C18, 3.0 x 30 mm, 2.7 µm, 600 bar		

## Checkout Samples

### Checkout Sample for UV Detection at 600 bar or higher

The 5188-6529 (RRLC checkout sample) serves as standard for systems with 600 bar or higher and contains 100 ng/μL each of nine components dissolved in water / acetonitrile (65/35). The nine components are:

- Acetanilide
- Acetophenone
- Propiophenone
- Butyrophenone
- Benzophenone
- Valerophenone
- Hexanophenone
- Heptanophenone
- Octanophenone

### Checkout Sample for UV Detection at Maximum 400 bar and FL Detection

The 01080-68704 (Isocratic standard) serves as standard for isocratic systems, systems with a FLD as only detector, and for systems with 400 bar pumps. It contains each of four components dissolved in methanol.

- Dimethylphthalate
- Diethylphthalate
- Biphenyl
- o-Terphenyl

## Checkout Sample for ELSD

The 5190-0488 (Caffeine standards kit) serves as standard for several instrument related procedures. The 200 µg/mL standard is used for checkout of ELSD systems. The standard kit contains calibrated amounts of caffeine in water with concentrations of:

- 0.5 µg/mL
- 1 µg/mL
- 2 µg/mL
- 5 µg/mL
- 25 µg/mL
- 50 µg/mL
- 100 µg/mL
- 200 µg/mL



## Checkout Method for Isocratic, 400 bar and FLD only systems

This checkout method covers configurations with isocratic or 400 bar pumps or systems with FLD as only detectors. For column options and checkout samples, see [Table 4 Overview of column options for different pumps](#) on page 70, and [Checkout Sample for UV Detection at Maximum 400 bar and FL Detection](#) on page 71.

### 1 Section NOT Applicable

2 Install the checkout column.

3 Setup the system.

a Set parameters for the pump.

OR

**Table 5:** Checkout method parameter settings G7111A/B, G7112B, G7120A, G7132A, G7104A/C, G7131A/C, or G5654A

Parameter	Value
Flow	1 mL/min (0.8 mL/min for G5654A)
Solvent A	Water
Solvent B	ACN
Compressibility	Use solvent types (Use 95 for G7111A/B) for Compressibility
Composition	35 % A (Water)
Composition	65 % B (ACN)
Stoptime	10 min
Pressure Limit	400 bar
Minimum Stroke	Automatic

- b** Set parameters for the injector.

**Table 6:** Checkout method parameter settings G7129A/B, G7167A/B, G7137A, or G5668A

Parameter	Value
Injection	1 µL
Stoptime	as pump
Draw speed	100 µL/min

OR

- c** Set parameters for the multicolumn thermostat.

**Table 7:** Checkout method parameter settings G7116A/B, or G7130A

Parameter	Value
Temperature (left)	40 °C
Temperature (right)	combined
Stoptime	as pump

- d** Set parameters for the detector.

**Table 8:** Checkout method parameter settings G7115A, G7165A, or G7117A/B/C

Parameter	Value
Signal A	254 /4 nm
Ref A	360 /100 nm
Peakwidth	40 Hz
Stoptime	as pump
Spectrum	None
Autobalance	Prerun

OR

**Table 9:** Checkout method parameter settings G7114A/B

Parameter	Value
Wavelength	254 nm
Peakwidth	40 Hz

**Quick Start Guide**

Installation Checkout (Customer presence recommended)

Parameter	Value
Stoptime	as pump
Autobalance	Prerun

OR

- 4 Start the system.
- 5 Equilibrate the system until the pressure signal and the detector baseline are stable.
- 6 Run the checkout sample and check the chromatogram for obvious abnormalities.

## Checkout Method for 1260 and 1290 Systems with UV Detectors

This checkout method covers configurations with gradient pumps with pressures of 600 bar and above with UV detectors. For column options and checkout samples, see [Table 4 Overview of column options for different pumps](#) on page 70 and [Checkout Sample for UV Detection at 600 bar or higher](#) on page 71.

### 1 Section NOT Applicable

2 Install the checkout column.

3 Setup the system.

a Set parameters for the pump.

**Table 10:** Checkout method parameter settings G7111A

Parameter	Value
Flow	1 mL/min
Solvent A	Water
Solvent B	ACN
Compressibility	75
Composition	35 % A (Water)
Composition	65 % B (ACN)
Stoptime	10 min
Pressure Limit	400 bar
Minimum Stroke	Automatic

OR

**Table 11:** Checkout method parameter settings G7111B, G7112B, G7120A, G7132A, G7104A/C, G7131A/C, or G5654A

Parameter	Value
Flow	0.8 mL/min (0.6 mL/min for G5654A)
Solvent A	Water
Solvent B	ACN
Compressibility	Use solvent types (use 95 for G7111B)

Parameter	Value
Composition	60 % A (Water)
Composition	40 % B (ACN)
Stoptime	10 min
Minimum Stroke	Automatic
Timetable	2.5 min, 80 %B

**b** Set parameters for the injector.

**Table 12:** Checkout method parameter settings G7129A/B, G7167A/B, G7137A, or G5668A

Parameter	Value
Injection	1 µL
Stoptime	as pump
Draw speed	100 µL/min

OR

**Table 13:** Checkout method parameter settings G7167C

Parameter	Value
Injection	1 µL
Stoptime	as pump
Draw speed	100 µL/min
<b>Feed Injection Mode</b>	<b>mandatory</b>
Feed Speed	Adaptive: 80 % of the pump flow
Flush-out	automatic
<b>Flow-through Injection Mode</b>	<b>optional</b> <sup>1</sup>
Flush-out Factor	3.5
Delay Volume Reduction	Enabled

<sup>1</sup> Flow-through Injection is optional. It can be done as an extra to the Feed Injection by customer request.

- c Set parameters for the multicolumn thermostat.

**Table 14:** Checkout method parameter settings G7116A/B, or G7130A

Parameter	Value
Temperature (left)	40 °C
Temperature (right)	combined
Stoptime	as pump

- d Set parameters for the multicolumn thermostat.

**Table 15:** Checkout method parameter settings G7116A/B, or G7130A

Parameter	Value
Temperature (left)	40 °C
Temperature (right)	combined
Stoptime	as pump

- e Set parameters for the detector.

**Table 16:** Checkout method parameter settings G7115A, G7165A, or G7117A/B/C

Parameter	Value
Signal A	254 /4 nm
Ref A	360 /100 nm
Peakwidth	40 Hz
Stoptime	as pump
Spectrum	None
Autobalance	Prerun

OR

**Table 17:** Checkout method parameter settings G7114A/B

Parameter	Value
Wavelength	254 nm
Peakwidth	40 Hz
Stoptime	as pump
Autobalance	Prerun

**Quick Start Guide**

Installation Checkout (Customer presence recommended)

- 4 Start the system.
- 5 Equilibrate the system until the pressure signal and the detector baseline are stable.
- 6 Run the checkout sample and check the chromatogram for obvious abnormalities.

## Checkout Method for 1260 and 1290 Systems with ELSD or RID

This checkout method covers configurations with RI or ELS detectors as only detectors. For column options and checkout samples (keep in mind the samples are different for ELSD and RID), see [Table 4 Overview of column options for different pumps](#) on page 70, [Checkout Sample for ELSD](#) on page 72, and [Checkout Sample for RID](#).

### 1 Section NOT Applicable

2 Install the restriction capillary in the MCT.

3 Setup the system.

a Set parameters for the pump.

**Table 18:** Checkout method parameter settings G7110B, G7111A/B, G7112B, G7120A, G7132A, G7104A/C, G7131A/C, or G5654A with ELSD or RID

Parameter	Value
Flow	1 mL/min
Solvents	Water
Compressibility	46
Stoptime	5 min
Pressure Limit	400 bar
Minimum Stroke	Automatic

b Set parameters for the injector.

**Table 19:** Checkout method parameter settings G7129A/B, G7167A/B, G7137A, or G5668A with ELSD or RID

Parameter	Value
Injection	20 µL
Stoptime	as pump
Draw speed	100 µL/min

OR



**Table 20:** Checkout method parameter settings G7167C

Parameter	Value
Injection	20 µL
Stoptime	as pump
Draw speed	100 µL/min
<b>Feed Injection Mode</b>	<b>mandatory</b>
Feed Speed	Adaptive: 80 % of the pump flow
Flush-out	automatic
<b>Flow-through Injection Mode</b>	<b>optional</b> <sup>1</sup>
Flush-out Factor	3.5
Delay Volume Reduction	Enabled

<sup>1</sup> Flow-through Injection is optional. It can be done as an extra to the Feed Injection by customer request.

**c** Set parameters for the multicolumn thermostat.

**Table 21:** Checkout method parameter settings G7116A/B with ELSD, RID

Parameter	Value
Temperature (left)	35 °C
Temperature (right)	combined
Stoptime	as pump

**d** Set parameters for the detector.

**Table 22:** Checkout method parameter settings G7102A, G426XA/B

Parameter	Value
Nebulizer	Temperature: 70 °C
Evaporator	Temperature: 70 °C
Gas Flow	1.60 SLM
Data Rate	10 Hz
Smoothing	1 s
PMT Gain	1 (No PMT Gain setting for G7102A)
LED Intensity	100 % (No Laser setting for the Agilent G7102A)

OR

**Table 23:** Checkout method parameter settings G7162A/B

Parameter	Value
Optical Unit Temperature	35 °C
Signal	Acquire
Response time	4 s
Signal Polarity	positive
Automatic Zero	on
Automatic Recycling	off

- 4 Start the system.
- 5 Equilibrate the system until the pressure signal and the detector baseline are stable.
- 6 Run the checkout sample and check the chromatogram for obvious abnormalities.



## 5 Parts and Consumables

This chapter provides information on additional parts and consumables.

### **HPLC System Tool Kit 84**

#### **InfinityLab Quick Connect and Quick Turn Fittings 85**

InfinityLab Quick Connect Fittings 85

InfinityLab Quick Connect Fitting Replacement Capillaries 87

InfinityLab Quick Turn Fitting 88

Capillaries for Use with the InfinityLab Quick Turn Fitting 89

#### **Safety Caps and Solvent Bottles 90**

Stay Safe Caps 90

Solvent Bottles 93

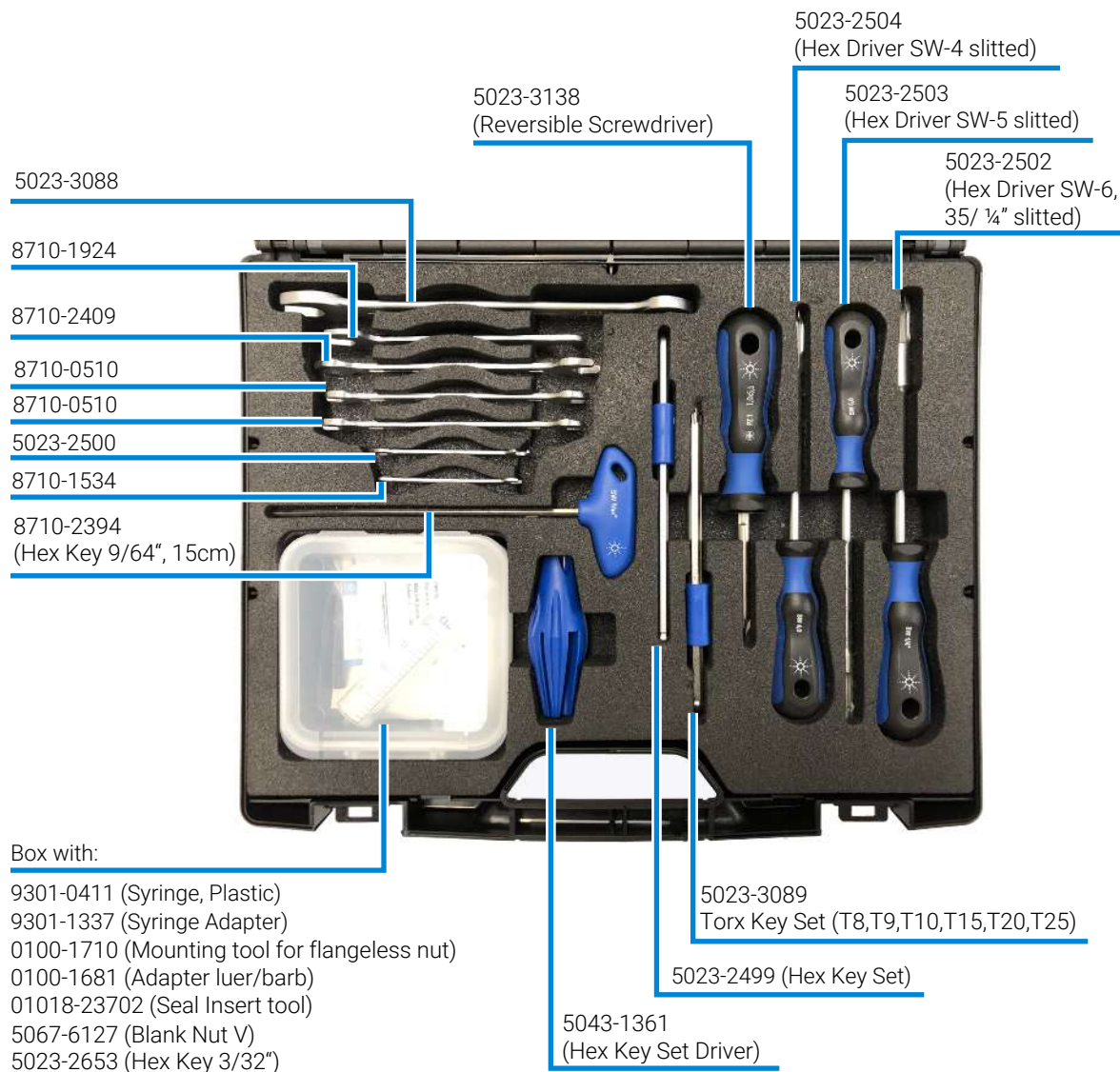
Additional Parts 94

#### **InfinityLab Flex Bench Family 95**

InfinityLab Flex Bench 95

InfinityLab Benchtop 96

## HPLC System Tool Kit



# InfinityLab Quick Connect and Quick Turn Fittings

## InfinityLab Quick Connect Fittings

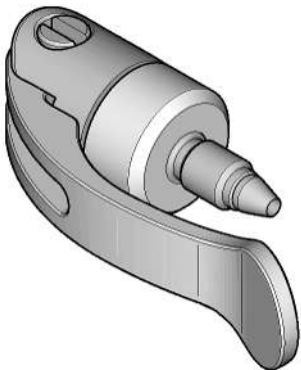


































Figure 24: InfinityLab Quick Connect Fitting

p/n	Description
 5067-5965	InfinityLab Quick Connect LC fitting (fitting without preinstalled capillary)
 5043-0924	Front Ferrule for Quick Connect/Turn Fitting
 5067-5961	InfinityLab Quick Connect Assy ST 0.075 mm x 105 mm
 5067-6163	InfinityLab Quick Connect Assy ST 0.075 mm x 150 mm
 5067-6164	InfinityLab Quick Connect Assy ST 0.075 mm x 220 mm
 5067-6165	InfinityLab Quick Connect Assy ST 0.075 mm x 280 mm
 5067-5957	InfinityLab Quick Connect Assy ST 0.12 mm x 105 mm
 5067-5958	InfinityLab Quick Connect Assy ST 0.12 mm x 150 mm
 5067-5959	InfinityLab Quick Connect Assy ST 0.12 mm x 220 mm
 5067-5960	InfinityLab Quick Connect Assy ST 0.12 mm x 280 mm
 5067-6166	InfinityLab Quick Connect Assy ST 0.17 mm x 105 mm
 5067-6167	InfinityLab Quick Connect Assy ST 0.17 mm x 150 mm

**Parts and Consumables**  
InfinityLab Quick Connect and Quick Turn Fittings

p/n	Description
 <a href="#">5067-6168</a>	InfinityLab Quick Connect Assy ST 0.17 mm x 220 mm
 <a href="#">5067-6169</a>	InfinityLab Quick Connect Assy ST 0.17 mm x 280 mm

## InfinityLab Quick Connect Fitting Replacement Capillaries

p/n	Description
 5500-1174	InfinityLab Capillary ST 0.075 mm x 105 mm
 5500-1175	InfinityLab Capillary ST 0.075 mm x 150 mm
 5500-1176	InfinityLab Capillary ST 0.075 mm x 220 mm
 5500-1177	InfinityLab Capillary ST 0.075 mm x 250 mm
 5500-1178	InfinityLab Capillary ST 0.075 mm x 280 mm
 5500-1173	InfinityLab Capillary ST 0.12 mm x 105 mm
 5500-1172	InfinityLab Capillary ST 0.12 mm x 150 mm
 5500-1171	InfinityLab Capillary ST 0.12 mm x 220 mm
 5500-1170	InfinityLab Capillary ST 0.12 mm x 280 mm
 5500-1179	InfinityLab Capillary ST 0.12 mm x 400 mm
 5500-1180	InfinityLab Capillary ST 0.12 mm x 500 mm
 5500-1181	InfinityLab Capillary ST 0.17 mm x 105 mm
 5500-1182	InfinityLab Capillary ST 0.17 mm x 150 mm
 5500-1183	InfinityLab Capillary ST 0.17 mm x 220 mm
 5500-1230	InfinityLab Capillary ST 0.17 mm x 280 mm
 5500-1231	InfinityLab Capillary ST 0.17 mm x 500 mm
 5500-1259	InfinityLab Capillary ST 0.25 mm x 150 mm
 5500-1260	InfinityLab Capillary ST 0.25 mm x 400 mm

InfinityLab Quick Turn Fitting

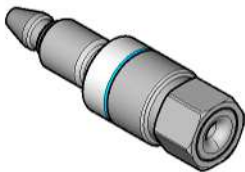




Figure 25: InfinityLab Quick Turn Fitting

p/n	Description
 <a href="#">5067-5966</a>	InfinityLab Quick Turn Fitting
 <a href="#">5043-0924</a>	Front Ferrule for Quick Connect/Turn Fitting



## Capillaries for Use with the InfinityLab Quick Turn Fitting

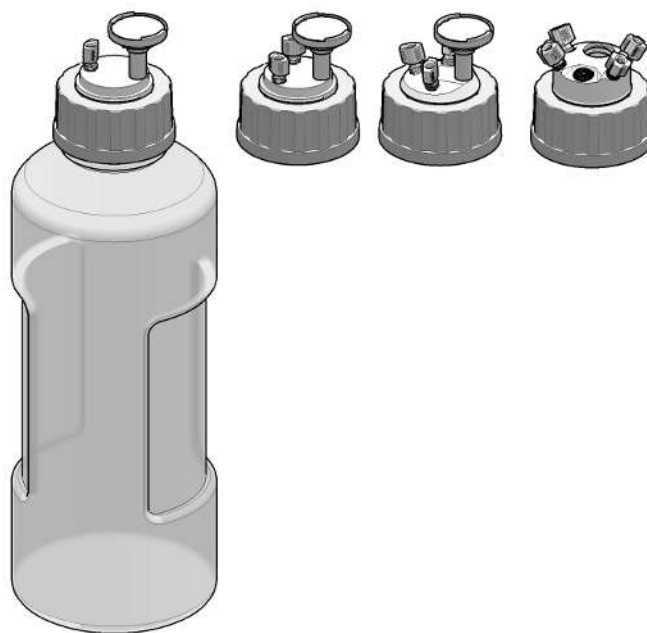
p/n	Description
 5500-1198	Capillary ST 0.075 mm x 105 mm, long socket
 5500-1232	Capillary ST 0.075 mm x 150 mm, long socket
 5500-1206	Capillary ST 0.075 mm x 250 mm, long socket
 5500-1205	Capillary ST 0.075 mm x 500 mm, long socket
 5500-1188	Quick Turn Capillary ST 0.12 mm x 105 mm, long socket
 5500-1189	Capillary ST 0.12 x 150 mm, long socket
 5500-1233	Capillary ST 0.12 mm x 180 mm, long socket
 5500-1190	Capillary ST 0.12 mm x 200 mm, long socket
 5500-1191	InfinityLab Quick Turn Capillary ST 0.12 mm x 280 mm, long socket
 5500-1192	Capillary ST 0.12 mm x 500 mm, long socket
 5500-1193	InfinityLab Quick Turn Capillary ST 0.17 mm x 105 mm, long socket
 5500-1194	Capillary ST 0.17 mm x 150 mm, long socket
 5500-1234	Capillary ST 0.17 mm x 180 mm
 5500-1195	Capillary ST 0.17 mm x 200 mm, long socket
 5500-1196	Capillary ST 0.17 mm x 280 mm, long socket
 5500-1235	Capillary ST 0.17 mm x 380 mm, long socket
 5500-1236	Capillary ST 0.17 mm x 400 mm, long socket
 5500-1197	Capillary ST 0.17 mm x 500 mm, long socket
 5500-1237	Capillary 0.17 mm x 700 mm, ns/ns
 5500-1262	Capillary 0.25 mm x 150 mm, ns/ns
 5500-1263	Capillary ST 0.25 mm x 400 mm, long socket
 5500-1200	Quick Turn Capillary ST 0.12 mm x 130 mm SL/M
 5500-1288	Capillary ST 0.12 mm x 150 mm, long socket, M4
 5500-1290	Capillary ST 0.17 mm x 150 mm, long socket, M4

## Safety Caps and Solvent Bottles

### Stay Safe Caps


**Table 24:** Stay Safe Caps

Part No.	Description	Fittings	Vent Ports	Filter Ports	Waste Ports
5043-1217	GL45 with 1 port 1 vent valve with time strip	1 x 3.2 mm	1		
5043-1218	GL45 with 2 ports 1 vent valve with time strip	2 x 3.2 mm	1		
5043-1219	GL45 with 3 ports 1 vent valve with time strip	3 x 3.2 mm	1		
5043-1220	GL45 with 4 ports 1 leak hose	4 (2 x 3.2 mm, 1 x 2.3 mm, 1 x 1.6 mm)		1	1



**Figure 26:** Solvent bottle with different types of Stay Safe caps

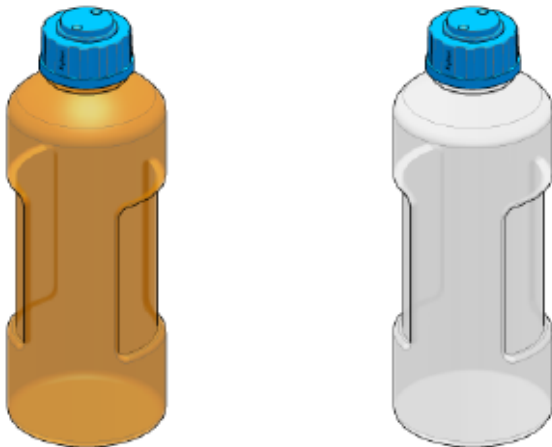
### Kits

p/n	Description
 <a href="#">5043-1221</a>	6 L waste can with 1 Stay Safe cap GL45 with 4 ports











**Figure 27:** Kit: Waste can with Stay Safe cap

## Solvent Bottles







**Figure 28:** Solvent bottles


p/n	Description
 9301-6523	Solvent bottle, clear, 500 mL with cap
 9301-6524	Solvent bottle, clear, 1000 mL
 9301-6525	Solvent bottle, amber, 500 mL with cap
 9301-6526	Solvent bottle, amber, 1000 mL
 9301-6527	Solvent bottle, clear, 125 mL
 9301-6528	Solvent bottle, clear, 1000 mL with cap
 9301-6529	Identification silicone ring (8/pk with 4 different colors)
 9301-6530	Sticker for solvent bottles (100/pk)

## Additional Parts

For fitting ports:

p/n	Description
 5043-1216	Fitting for 3.2 mm tubing, PFA, 2/pk
 5043-1215	Fitting for 2.3 mm tubing, PFA, 2/pk
 5043-1214	Fitting for 1.6 mm tubing, PFA, 2/pk
 5043-1198	Screw plug 1/8 in, PTFE, 2/pk


For vent port:

p/n	Description
 5043-1190	Venting valve with time strip, PTFE, 1 µm



For filter port:

p/n	Description
 5043-1193	Charcoal filter with time strip (58 g) for waste container

For waste port:



p/n	Description
 5043-1207	2-ports waste collector, PTFE

Miscellaneous:

p/n	Description
 5043-1191	Thread adapter PTFE GL45 (M) - GL40 (F)
 5043-1192	Thread adapter PTFE GL45 (M) - GPI 38-430 (F)






# InfinityLab Flex Bench Family

## InfinityLab Flex Bench



p/n	Description
 5043-1252	InfinityLab Flex Bench
 5043-1759	InfinityLab Flex Bench with pre-assembled power strip

The Flex Bench includes the framework, four shelf assemblies, and one waste bin.

Accessories:





p/n	Description
 5043-1287	Shelf assembly
 8121-1245	Valve shelf assembly
 5043-1278	Waste bin
 8121-2258	Power cord
 5043-1289	InfinityLab Flex Bench Replacement Kit (spare parts: screws and casters)

InfinityLab Benchtop

p/n	Description
 <a href="#">5043-1711</a>	InfinityLab Benchtop
 <a href="#">5043-1740</a>	InfinityLab Benchtop with pre-assembled power strip

The Benchtop includes the framework and three shelf assemblies.

Accessories:

p/n	Description
 <a href="#">5043-1750</a>	Shelf assembly
 <a href="#">8121-1245</a>	Valve shelf assembly
 <a href="#">8121-2258</a>	Power cord
 <a href="#">5043-1289</a>	InfinityLab Flex Bench Replacement Kit (spare parts: screws and casters)



## In This Book

This manual contains technical reference information about the Agilent 1290 Infinity III LC System.

The manual describes the following:

- introduction,
- install the modules,
- configuration settings,
- quick start guide,
- parts and consumables.

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies Inc. 2014-2025  
Edition: 05/2025

Document No: SD-29000160 Rev. E.01

