



Agilent 1260 Infinity Bio-inert Analytical-scale Fraction Collector

User Manual



BIO
inert



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2011-2012

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.

Manual Part Number

G5664-90001 Rev. B

Edition

05/2012

Printed in Germany

Agilent Technologies
Hewlett-Packard-Strasse 8
76337 Waldbronn

This product may be used as a component of an in vitro diagnostic system if the system is registered with the appropriate authorities and complies with the relevant regulations. Otherwise, it is intended only for general laboratory use.

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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will

receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

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.

In This Guide

This manual contains technical reference information about the Agilent 1260 Infinity Bio-inert Analytical-scale Fraction Collector G5664A.

1 Introduction

This chapter gives an introduction to the module..

2 Site Requirements and Specifications

This chapter provides information on environmental requirements, physical and performance specifications.

3 Installing the Fraction Collector

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

4 Using the Fraction Collector

This chapter provides information on how to set up the module.

5 How to Optimize the Fraction Collector

This chapter provides information on how to optimize the module.

6 Troubleshooting and Diagnostics

Overview about the troubleshooting and diagnostic features.

7 Error Information

This chapter describes the meaning of error messages, and provides information on probable causes and suggested actions how to recover from error conditions.

8 Test Functions

This chapter describes the module's built in test functions.

9 Maintenance

This chapter describes the maintenance of the module.

10 Parts for Maintenance

This chapter provides information on parts for maintenance.

11 Identifying Cables

This chapter provides information on cables used with the Agilent modules.

12 Hardware Information

This chapter describes the module in more detail on hardware and electronics.

13 Appendix

This chapter provides addition information on safety, legal and web.

Contents

1	Introduction	9
	Introduction to the Fraction Collector	10
	Bio-inert Materials	10
	Early Maintenance Feedback	12
	Instrument Layout	13
2	Site Requirements and Specifications	15
	Site Requirements	16
	Physical Specifications	19
	Performance Specifications	20
3	Installing the Fraction Collector	23
	Unpacking the module	24
	Optimizing the Stack Configuration	26
	Installing the Module	30
	Installing a Thermostatted Fraction Collector	35
	Flow Connections to the Fraction Collector	40
	Fraction Collector Trays	43
	Configure Well-plate Types	46
	Transporting the Fraction Collector	50
4	Using the Fraction Collector	51
	Configuration of the Fraction Collector	52
	Delay Volumes and Delay Calibration	58
	Setting up a Fraction Collector Method	63
	Starting Your Run with Fraction Collection	69
	Viewing Your Results	74
	Special Applications	77
	Limitations and How to Avoid Problems	80
	Check-out Procedures	81
	Application Notes	84
	Solvent Information	85

5	How to Optimize the Fraction Collector	89
	Optimizing Fraction Collection	90
6	Troubleshooting and Diagnostics	91
	Overview of the Module's Indicators and Test Functions	92
	Status Indicators	93
	Agilent Lab Advisor Software	95
7	Error Information	97
	What Are Error Messages	99
	General Error Messages	100
	Module Error Messages	106
8	Test Functions	117
	Maintenance Functions	118
	Transport Unit Self Alignment	120
	Step Commands	122
9	Maintenance	125
	Introduction to Maintenance	126
	Cautions and Warnings	127
	Cleaning the Module	129
	Overview of Maintenance	130
	Replacing the Inlet/Waste Tubings	132
	Replacing the Valve to Needle Tubing	136
	Exchanging the Needle Assembly	140
	Exchanging the Needle/Capillary Carrier Assembly	143
	Exchanging the Diverter Valve	148
	Exchanging the Internal Tray	152
	Repairing or Exchanging a Funnel of the Internal Tray	154
	Exchanging the Leak Sensor	157
	Exchanging or Installing the (Optional) Interface Board	160
	Replacing Module Firmware	162

10 Parts for Maintenance 163

Fraction Collector Main Assemblies	164
Supported Trays for a Fraction Collector	166
List of Recommended Vials and Caps	167
Recommended Plates and Closing Mats	170
Fraction Collector Accessory Kit	172
Fraction Collector Thermostat	173
Needle Assembly	174
Diverter-Valve Assembly	175
Tubing Kits	176
Internal Tray Assembly	177

11 Identifying Cables 179

Cable Overview	180
Analog Cables	182
Remote Cables	184
BCD Cables	187
CAN/LAN Cables	189
External Contact Cable	190
RS-232 Cable Kit	191

12 Hardware Information 193

Electrical Connections	194
Interfaces	196
Setting the 8-bit Configuration Switch (without On-board) LAN	202
Optional Interface Boards	206

13 Appendix 211

General Safety Information	212
The Waste Electrical and Electronic Equipment Directive	215
Lithium Batteries Information	216
Radio Interference	217
Sound Emission	218
Agilent Technologies on Internet	219



1

Introduction

Introduction to the Fraction Collector 10

Bio-inert Materials 10

Early Maintenance Feedback 12

Instrument Layout 13

This chapter gives an introduction to the module.



Introduction to the Fraction Collector

The G5664A is a fraction collector for analytical scale fractionation. The bio-inert fraction collector is designed with a metal free flow path for the sample. The pH range is from 1 – 13 (short term 14 pH).

Bio-inert Materials

For the Agilent 1260 Infinity Bio-inert LC system, Agilent Technologies uses highest quality materials in the flow path (also referred to as wetted parts), which are widely accepted by life scientists, as they are known for optimum inertness to biological samples and ensure best compatibility with common samples and solvents over a wide pH range. Explicitly, the complete flow path is free of stainless steel and free of other alloys containing metals such as iron, nickel, cobalt, chromium, molybdenum or copper, which can interfere with biological samples. The flow downstream of the sample introduction contains no metals whatsoever.

Table 1 Bio-inert Materials

Module	Materials
Agilent 1260 Infinity Bio-inert Quaternary Pump (G5611A)	Titanium, gold, platinum-iridium, ceramic, ruby, PTFE, PEEK
Agilent 1260 Infinity Bio-inert High-Performance Autosampler (G5667A)	Upstream of sample introduction: • Titanium, gold, PTFE, PEEK, ceramic Downstream of sample introduction: • PEEK, ceramic
Agilent 1260 Infinity Bio-inert Manual Injector (G5628A)	PEEK, ceramic
Agilent 1260 Infinity Bio-inert Analytical Fraction Collector (G5664A)	PEEK, ceramic, PTFE

Table 1 Bio-inert Materials

Module	Materials
Bio-inert Flow Cells:	
Standard flow cell bio-inert, 10 mm, 13 µl, 120 bar (12 MPa) for MWD/DAD, includes Capillary Kit Flow Cells BIO (p/n G5615-68755) (p/n G5615-60022) (for Agilent 1260 Infinity Diode Array Detectors DAD G1315C/D)	PEEK, ceramic, sapphire, PTFE
Max-Light Cartridge Cell Bio-inert (10 mm, V(σ) 1.0 µl) (p/n G5615-60018) and Max-Light Cartridge Cell Bio-inert (60 mm, V(σ) 4.0 µl) (p/n G5615-60017) (for Agilent 1200 Infinity Series Diode Array Detectors DAD G4212A/B)	PEEK, fused silica
Bio-inert flow cell, 8 µL, 20 bar (pH 1–12) includes Capillary Kit Flow Cells BIO (p/n G5615-68755) (p/n G5615-60005) (for Agilent 1260 Infinity Fluorescence Detector FLD G1321B)	PEEK, fused silica, PTFE
Bio-inert heat-exchanger G5616-60050 for 1290 Infinity Thermostatted Column Compartment (G1316C)	PEEK (steel-cladded)
Bio-inert Valve heads	G4235A, G5631A, G5639A: PEEK, ceramic (Al ₂ O ₃ based)
Bio-inert Connection capillaries	Upstream of sample introduction: • Titanium Downstream of sample introduction: • Agilent uses stainless-steel-cladded PEEK capillaries, which keep the flow path free of steel and provide pressure stability to more than 600 bar.

NOTE

To ensure optimum bio-compatibility of your Agilent 1260 Infinity Bio-inert LC system, do not include non-inert standard modules or parts to the flow path. Do not use any parts that are not labeled as Agilent “Bio-inert”. For solvent compatibility of these materials, see [“Solvent information for parts of the 1260 Infinity Bio-inert LC system”](#) on page 85.

Early Maintenance Feedback

Maintenance requires the exchange of components which are subject to wear or stress. Ideally, the frequency at which components are exchanged should be based on the intensity of usage of the module and the analytical conditions, and not on a predefined time interval. The early maintenance feedback (**EMF**) feature monitors the usage of specific components in the instrument, and provides feedback when the user-selectable limits have been exceeded. The visual feedback in the user interface provides an indication that maintenance procedures should be scheduled.

EMF Counters

EMF counters increment with use and can be assigned a maximum limit which provides visual feedback in the user interface when the limit is exceeded. Some counters can be reset to zero after the required maintenance procedure.

Using the EMF Counters

The user-settable **EMF** limits for the **EMF Counters** enable the early maintenance feedback to be adapted to specific user requirements. The useful maintenance cycle is dependent on the requirements for use. Therefore, the definition of the maximum limits need to be determined based on the specific operating conditions of the instrument.

Setting the EMF Limits

The setting of the **EMF** limits must be optimized over one or two maintenance cycles. Initially the default **EMF** limits should be set. When instrument performance indicates maintenance is necessary, take note of the values displayed by the **EMF counters**. Enter these values (or values slightly less than the displayed values) as **EMF** limits, and then reset the **EMF counters** to zero. The next time the **EMF counters** exceed the new **EMF** limits, the **EMF** flag will be displayed, providing a reminder that maintenance needs to be scheduled.

Instrument Layout

The industrial design of the module incorporates several innovative features. It uses Agilent's E-PAC concept for the packaging of electronics and mechanical assemblies. This concept is based upon the use of expanded polypropylene (EPP) layers of foam plastic spacers in which the mechanical and electronic boards components of the module are placed. This pack is then housed in a metal inner cabinet which is enclosed by a plastic external cabinet. The advantages of this packaging technology are:

- virtual elimination of fixing screws, bolts or ties, reducing the number of components and increasing the speed of assembly/disassembly,
- the plastic layers have air channels molded into them so that cooling air can be guided exactly to the required locations,
- the plastic layers help cushion the electronic and mechanical parts from physical shock, and
- the metal inner cabinet shields the internal electronics from electromagnetic interference and also helps to reduce or eliminate radio frequency emissions from the instrument itself.

1 Introduction

Instrument Layout



2

Site Requirements and Specifications

Site Requirements 16

Physical Specifications 19

Performance Specifications 20

This chapter provides information on environmental requirements, physical and performance specifications.



Site Requirements

Power Considerations

The module power supply has wide ranging capability. It accepts any line voltage in the range described in [Table 2](#) on page 19. Consequently there is no voltage selector in the rear of the module. There are also no externally accessible fuses, because automatic electronic fuses are implemented in the power supply.

WARNING

Hazard of electrical shock or damage of your instrumentation can result, if the devices are connected to a line voltage higher than specified.

→ Connect your instrument to the specified line voltage only.

WARNING

The module is partially energized when switched off, as long as the power cord is plugged in.

Repair work at the module can lead to personal injuries, e.g. electrical shock, when the cover is opened and the module is connected to power.

→ Always unplug the power cable before opening the cover.

→ Do not connect the power cable to the instrument while the covers are removed.

CAUTION

Inaccessible power plug.

In case of emergency it must be possible to disconnect the instrument from the power line at any time.

→ Make sure the power connector of the instrument can be easily reached and unplugged.

→ Provide sufficient space behind the power socket of the instrument to unplug the cable.

Power Cords

Different power cords are offered as options with the module. The female end of all power cords is identical. It plugs into the power-input socket at the rear. The male end of each power cord is different and designed to match the wall socket of a particular country or region.

WARNING

Absence of ground connection or use of unspecified power cord

The absence of ground connection or the use of unspecified power cord can lead to electric shock or short circuit.

- Never operate your instrumentation from a power outlet that has no ground connection.
 - Never use a power cord other than the Agilent Technologies power cord designed for your region.
-

WARNING

Use of unsupplied cables

Using cables not supplied by Agilent Technologies can lead to damage of the electronic components or personal injury.

- Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.
-

WARNING

Unintended use of supplied power cords

Using power cords for unintended purposes can lead to personal injury or damage of electronic equipment.

- Never use the power cords that Agilent Technologies supplies with this instrument for any other equipment.
-

Bench Space

The module dimensions and weight (see [Table 2](#) on page 19) allow you to place the module on almost any desk or laboratory bench. It needs an additional 2.5 cm (1.0 inches) of space on either side and approximately 8 cm (3.1 inches) in the rear for air circulation and electric connections.

If the bench shall carry a complete HPLC system, make sure that the bench is designed to bear the weight of all modules.

The module should be operated in a horizontal position.

Condensation

CAUTION

Condensation within the module

Condensation will damage the system electronics.

- Do not store, ship or use your module under conditions where temperature fluctuations could cause condensation within the module.
 - If your module was shipped in cold weather, leave it in its box and allow it to warm slowly to room temperature to avoid condensation.
-

Physical Specifications

Table 2 Physical Specifications

Type	Specification	Comments
Weight	13.5 kg (29.8 lbs)	
Dimensions (height × width × depth)	200 x 345 × 440 mm (8 x 13.5 × 17 inches)	
Line voltage	100 – 240 VAC, ± 10 %	Wide-ranging capability
Line frequency	50 or 60 Hz, ± 5 %	
Power consumption	180 VA / 180 W / 239 BTU	Maximum
Ambient operating temperature	4 - 40 °C (41 - 104 °F)	
Ambient non-operating temperature	-40 – 70 °C (-4 – 158 °F)	
Humidity	< 95 %, at 25 – 40 °C (77 – 104 °F)	Non-condensing
Operating altitude	Up to 2000 m (6562 ft)	
Non-operating altitude	Up to 4600 m (15091 ft)	For storing the module
Safety standards: IEC, CSA, UL	Installation category II, Pollution degree 2	For indoor use only.

Performance Specifications

Table 3 Specifications - Agilent 1260 Infinity Analytical Bio-inert Fraction Collector (G5664A)

Type	Specifications
Delay volume	Approximately 50 μ L
Maximum system flow	10 mL/min
Fraction containers	Shallow or deep well plates up to 48 mm height (96 or 384 format) Test tubes up to 48 mm height Autosampler Vials (2 mL and 6 mL) Eppendorf safe-lock tubes (0.5 mL, 1.5 mL, 2.5 mL)
Fraction trays	Full trays (cover complete fraction collector) 4 well plates 40 x 20 mL test tubes (30 mm od, 48 mm height) 60 x 15 mL test tubes (25 mm od, 48 mm height) 126 x 8 mL test tubes (16 mm od, 48 mm height) 215 x 5 mL test tubes (12 mm od, 48 mm height) Trays with possibiity to add 1 additional half tray 2 well plates + 10 x 2 mL vials (+ one possible additional half tray) 100 x 2 mL vials (+ one possible additional half tray) Half trays (up to 3 per fraction collector) 15 x 6 mL vial 40 x 2 mL vial Plates for well plate trays (2 or 4 per fraction collector depending on well plate tray) Eppendorf safe-lock tubes (27 x 0.5 mL, 1.5 mL or 2.5 mL) 24 test tubes (18 mm OD) 54 x 2 mL vial 15 x 6 mL vial
Cooling	Optional
Trigger modes	Time slices and peak (threshold, up-/downslope, upper threshold and timetable), Boolean logic for different detector signals, combination of different modes, manual trigger (supported with Agilent 1200 Series Instant Pilot)

Table 3 Specifications - Agilent 1260 Infinity Analytical Bio-inert Fraction Collector (G5664A)

Type	Specifications
Trigger source	Agilent 1200 Infinity Series VWD, MWD and DAD detectors, Agilent 6100 Series Quadrupole LC/MS, ELSD, FLD, RID, third party detectors (require UIB)
Diverter valve	3/2 valve, switching time < 100 ms
Max. pressure	6 bar (diverter valve)
Environment	4 – 55 °C constant temperature, < 95 % humidity (non-condensing)
Materials in flow path	PEEK, Ceramic, PTFE
pH range	1 – 13 (short term 14) ¹

¹ For solvent compatibility, refer to section "Solvent information for parts of the 1260 Infinity Bio-inert LC system" in the manual

NOTE

Only one type of well-plates can be used at a time in one tray.

2 Site Requirements and Specifications

Performance Specifications



3

Installing the Fraction Collector

Unpacking the module	24
Optimizing the Stack Configuration	26
Installing the Module	30
Installing a Thermostatted Fraction Collector	35
Flow Connections to the Fraction Collector	40
Fraction Collector Trays	43
Configure Well-plate Types	46
Transporting the Fraction Collector	50

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



Unpacking the module

Damaged Packaging

If the delivery packaging shows signs of external damage, please call your Agilent Technologies sales and service office immediately. Inform your service representative that the instrument may have been damaged during shipment.

CAUTION

"Defective on arrival" problems

If there are signs of damage, please do not attempt to install the module. Inspection by Agilent is required to evaluate if the instrument is in good condition or damaged.

- Notify your Agilent sales and service office about the damage.
 - An Agilent service representative will inspect the instrument at your site and initiate appropriate actions.
-

Delivery Checklist

Ensure all parts and materials have been delivered with the fraction collector. For this compare the shipment content with the checklist included in each instrument box. Please report missing or damaged parts to your local Agilent Technologies sales and service office.

The following models of Agilent 1260 Infinity Bio-inert Fraction Collector are available:

- G5664A Bio-inert Fraction Collector, designed for flow rates below 10 mL/min, well-plates and a 40-funnel tray connecting to external locations of any size.
- The G5664A can also be equipped with a thermostat (G1330B).

Table 4 Delivery Checklist for the G5664A Bio-inert Analytical Scale Fraction Collector

Description	Quantity
Fraction collector module with inlet / waste tubing assembly	1
Power cord, local (Matching the sockets in your country or region)	1
Accessory kit, see “Fraction Collector Accessory Kit” on page 172	1
Delay sensor calibrant (p/n G1946-85020)	1

Table 5 Delivery Checklist for the G1330B Thermostat Module (optional)

Description	Quantity
Thermostat Module	1
Power cord, local (Matching the sockets in your country or region)	1
Accessory kit thermostat module (see “Fraction Collector Thermostat” on page 173)	1

Optimizing the Stack Configuration

If your fraction collector is part of a HPLC system, you can ensure optimum performance and minimum delay volume by installing the following configuration.

[Figure 1](#) on page 27 and [Figure 2](#) on page 28 show the configuration recommended for the fraction collector in an analytical scale system.

NOTE

Agilent 1260 Infinity LC Purification Systems can comprise many different LC modules. This section describes some common stack configurations and general considerations. On more complex systems, e.g. with two or more fraction collectors or with mass selective detector, other stack configurations might show advantages.

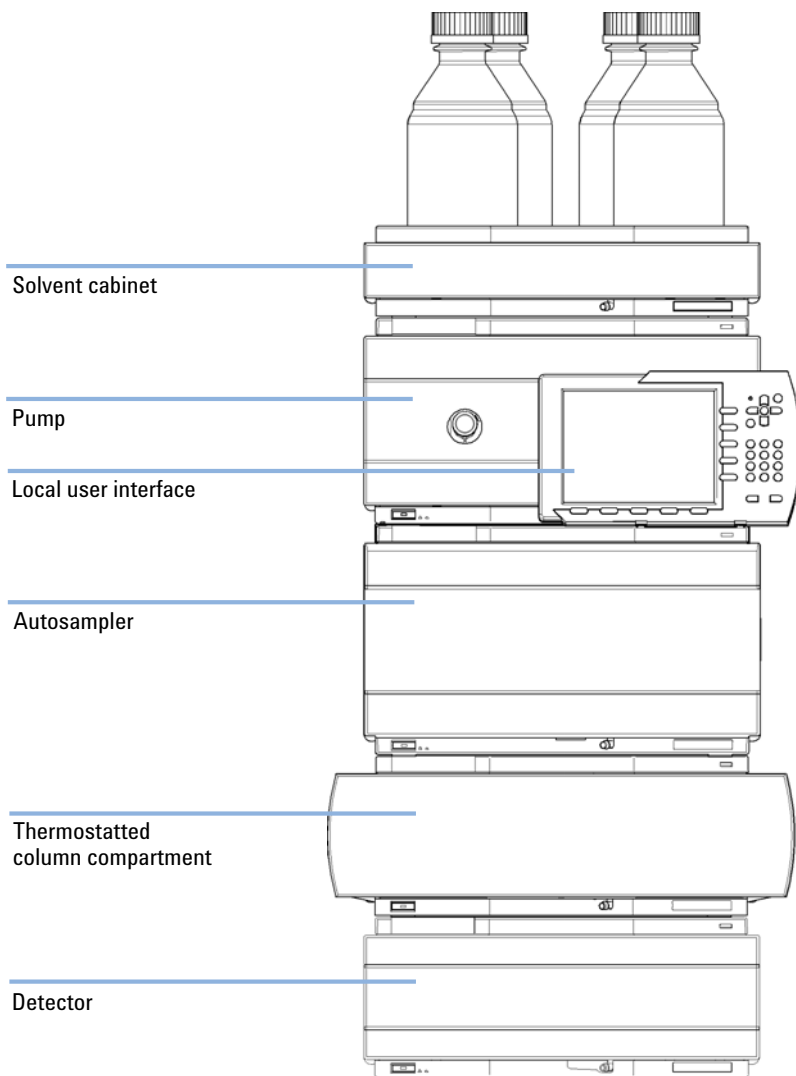


Figure 1 Recommended Stack Configuration (Front View)

3 Installing the Fraction Collector

Optimizing the Stack Configuration

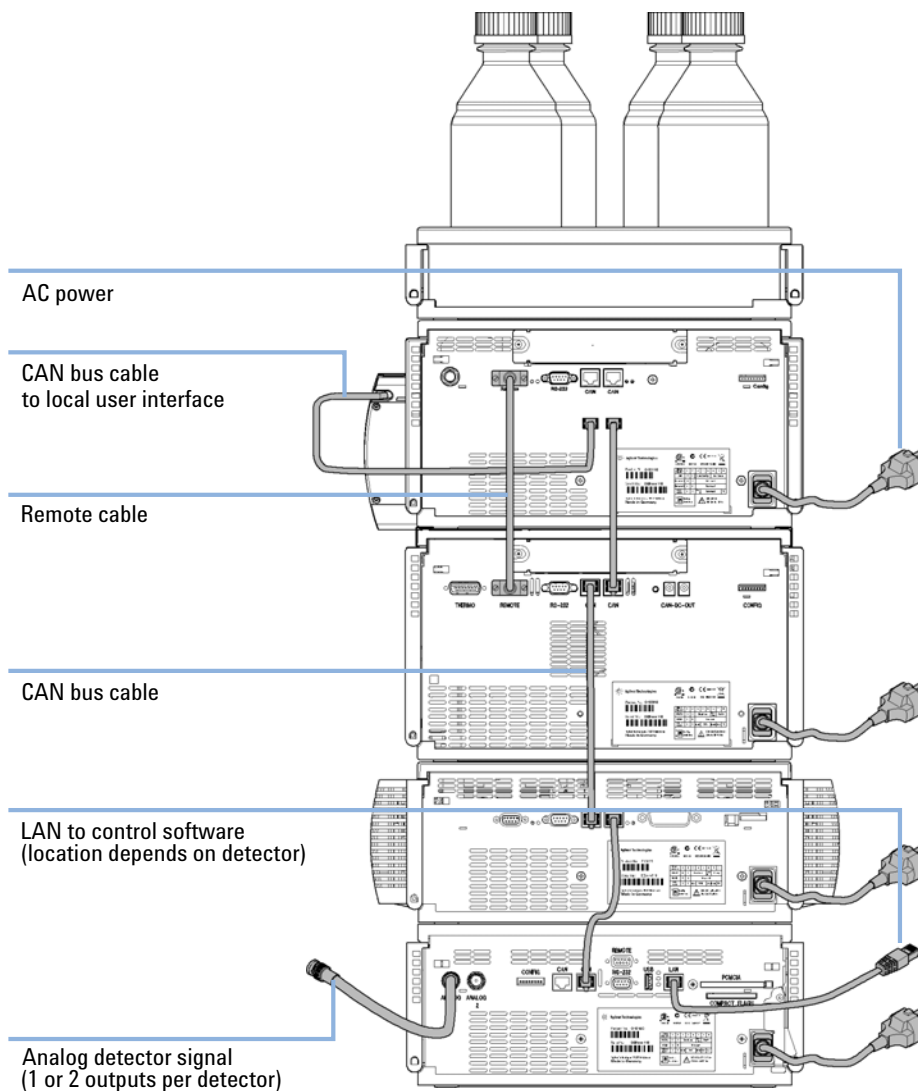


Figure 2 Recommended Stack Configuration (Rear View)

NOTE

If a G1330B ALS thermostat is part of the system it must always be installed underneath the thermostatted injector or fraction collector (see [“Installing a Thermostatted Fraction Collector”](#) on page 35).

NOTE

The fraction collector should never be positioned on top of a module that generates heat. This could lead to an unwanted evaporation of fractions in the fraction collector (e.g. Agilent 1260 Infinity Thermostatted Column Compartment G1316A or Agilent 1260 Infinity Diode Array Detectors).

Installing the Module

Parts required

Description

Fraction Collector
Chemstation or Agilent Instant Pilot G4208A
Power cord

For other cables see below and [“Cable Overview”](#) on page 180.

Preparations

Locate bench space
Provide power connections
Unpack the fraction collector

WARNING

When opening capillary or tube fittings solvents may leak out.

The handling of toxic and hazardous solvents and reagents can hold health risks.

→ Please observe appropriate safety procedures (for example, goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet supplied by the solvent vendor, especially when toxic or hazardous solvents are used.

WARNING

Personal injury

In the needle area there is a risk of being hurt by the needle during operation.

→ Keep fingers away from the needle area during fraction collector operation. Do not attempt to insert or remove a vial or a plate when the needle is positioned.

WARNING

The module is partially energized when switched off, as long as the power cord is plugged in.

Repair work at the module can lead to personal injuries, e.g. electrical shock, when the cover is opened and the module is connected to power.

- Always unplug the power cable before opening the cover.
 - Do not connect the power cable to the instrument while the covers are removed.
-

- 1 Install the LAN interface board in the fraction collector (if required), see [“Exchanging or Installing the \(Optional\) Interface Board”](#) on page 160.
- 2 Remove the adhesive tape which covers the side and front doors.
- 3 Open the front door and remove the left side door.
- 4 Remove the transport protection foam.
- 5 Install the corrugated waste tube in the plastic port at the front bottom center of the instrument. Slide the waste tubing coming from the internal tray (if present) through the plastic port and the corrugated waste tube (see

3 Installing the Fraction Collector

Installing the Module

Figure 3 on page 32). Route the corrugated waste tubing into a waste container.

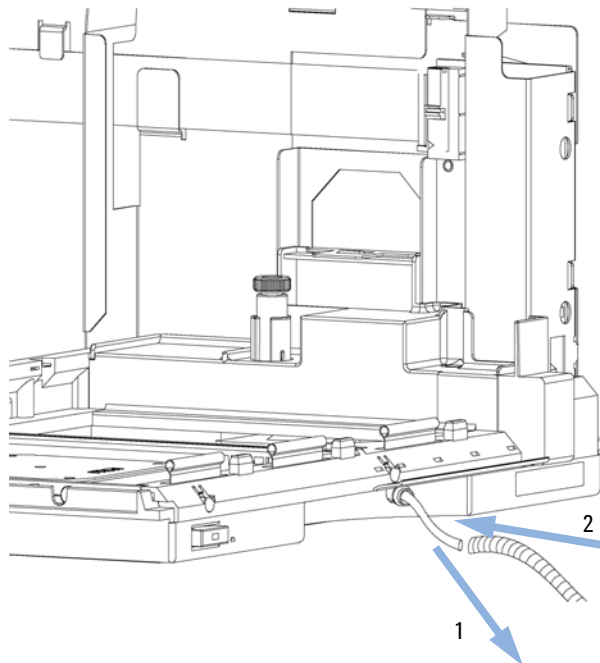


Figure 3 Installing the Corrugated Waste Tubing in the Plastic Port

- 6 Re-install the left side door (take care of the magnet at the back). Ensure the side door is correctly installed (its presence is sensed by a hall sensor, a missing side door will result in a NOT-READY state of the instrument).
- 7 Place the fraction collector in the stack or on the bench in all horizontal position.
- 8 Ensure the power switch at the front of the fraction collector is OFF.
- 9 Connect the power cable to the power connector at the rear of the fraction collector.

- 10** Connect the CAN cable to the other Agilent 1260 Infinity modules.
- 11** If an Agilent ChemStation is the controller, connect the LAN connection to the LAN interface (should be installed to the detector).
- 12** Connect the APG remote cable (optional) for non Agilent 1260 Infinity instruments.
- 13** Install the tray that has been delivered for your fraction collector. The test tube trays can be adjusted in height, depending on the height of the used test tubes. To adjust the height of the trays, press against the snappers at the guides in the 4 corners of a tray and move the top plate of the tray up or down (see [Figure 4](#) on page 33).

NOTE

The Agilent 1260 Infinity Bio-inert Analytical Scale Fraction Collector (G5664A) is shipped with two trays (40 x 2 ml half tray and 15 x 6 ml half tray). All other trays have to be ordered separately.

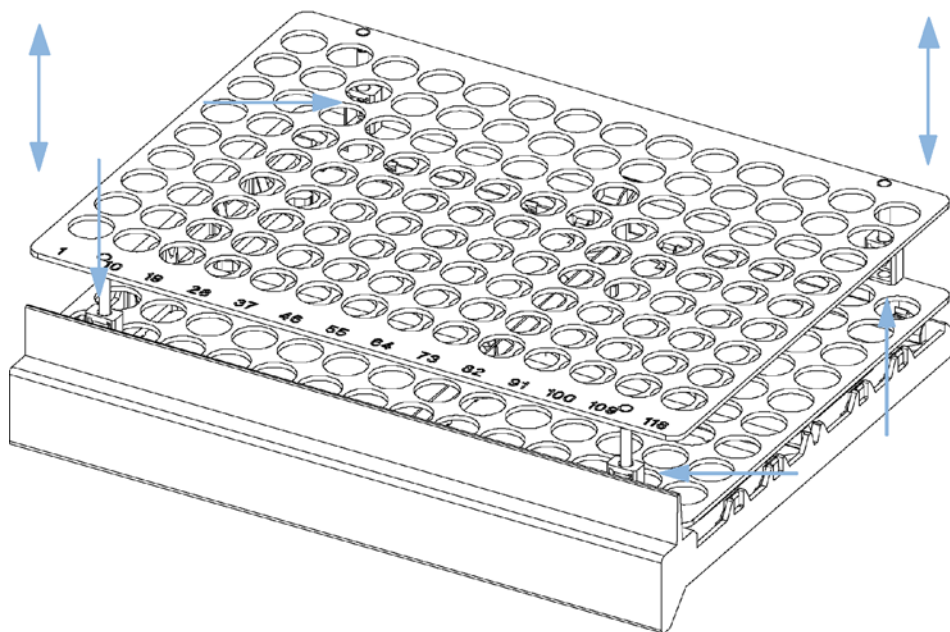


Figure 4 Adjusting the Height of the (Full) Test Tube

3 Installing the Fraction Collector

Installing the Module

- 14 Turn ON power by pushing the button at the lower left hand side of the fraction collector.
- 15 The exhaust fan will turn ON and remove potential solvent vapor from the inside of the instrument. After 2 minutes close the front door. Then the fraction collector will start the hardware initialization process. At the end of this process the status LED should be green.

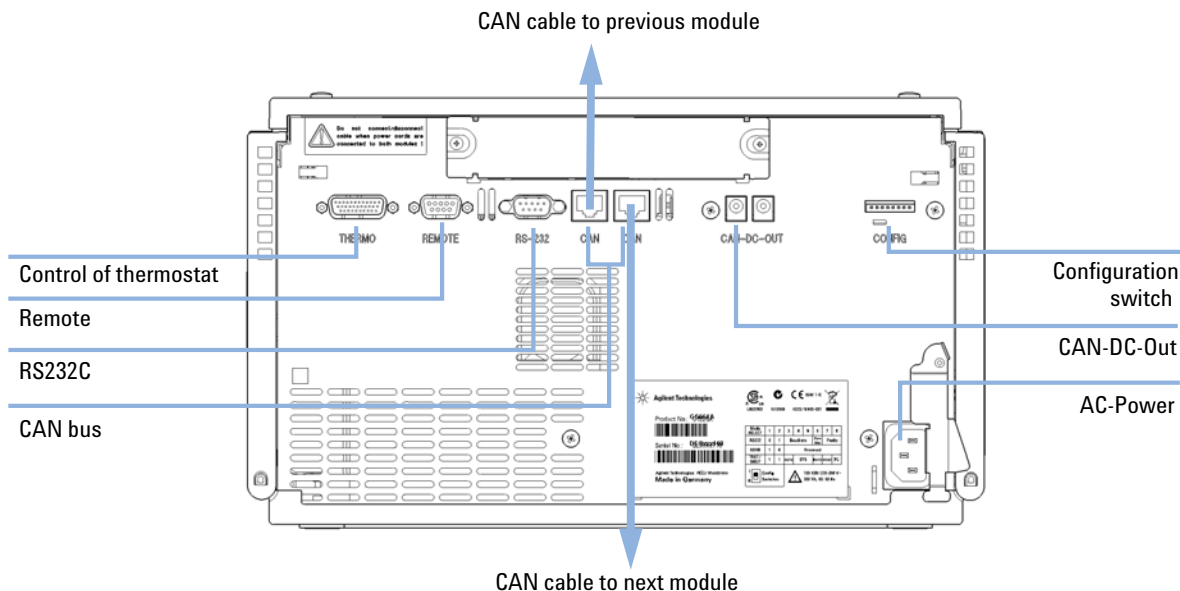


Figure 5 Cable Connections

NOTE

The fraction collector is turned ON when the line power switch is pressed and the green indicator lamp is illuminated. The detector is turned OFF when the line power switch is protruding and the green light is OFF.

Installing a Thermostatted Fraction Collector

Parts required	Description
	Fraction Collector and thermostat
	Chemstation or Agilent Instant Pilot G4208A
	Power cord

For other cables see below and [“Cable Overview”](#) on page 180.

Preparations	Locate bench space
	Provide power connections
	Unpack the fraction collector and the thermostat

WARNING

Personal injury

In the needle area there is a risk of being hurt by the needle during operation.

- Keep fingers away from the needle area during fraction collector operation. Do not attempt to insert or remove a vial or a plate when the needle is positioned.

WARNING

The module is partially energized when switched off, as long as the power cord is plugged in.

Repair work at the module can lead to personal injuries, e.g. electrical shock, when the cover is opened and the module is connected to power.

- Always unplug the power cable before opening the cover.
- Do not connect the power cable to the instrument while the covers are removed.

CAUTION

Damage of module electronics

Connecting or disconnecting the fraction collector to thermostat cable as long as the power cords are connected will damage the electronics of the modules.

- Before connecting the cable between the fraction collector and the thermostat, unplug the power cords of the fraction collector and the thermostat.

3 Installing the Fraction Collector

Installing a Thermostatted Fraction Collector

CAUTION

Damage through condensation

If the condensation tube is located in liquid the condensed water cannot flow out of the tube and the outlet is blocked. Any further condensation will then remain in the instrument. This may damage the instruments electronics.

→ Make sure that the condensation tube is always above the liquid level in the vessel.

CAUTION

Overheating of the module

Insufficient air circulation inside the module can lead to overheating.

→ Place the fraction collector thermostat with 25 cm (10 inch) space on each side for sufficient air circulation.

- 1 Place the thermostat on the bench.
- 2 Remove the front cover and route the condensation drain tube to a waste container.
- 3 Install the LAN interface board in the fraction collector (if required), see [“Exchanging or Installing the \(Optional\) Interface Board”](#) on page 160.
- 4 Remove the adhesive tape which covers the side and front doors.
- 5 Open the front door and remove the left side door.
- 6 Remove the transport protection foam.
- 7 Install the corrugated waste tube in the plastic port at the front bottom center of the fraction collector and route down into a waste container. Slide the waste tubing coming from the internal tray (if present) through the plastic port and the corrugated waste tube (see [Figure 3](#) on page 32). Route the corrugated waste tubing into a waste container.
- 8 Re-install the left side door (take care of the magnet at the back). Ensure the side door is correctly installed (its presence is sensed by a hall sensor, a missing side door will result in a NOT-READY state of the instrument).
- 9 Place the fraction collector on top of the thermostat. Make sure that the fraction collector is correctly engaged in the thermostat locks.

- 10** Remove the plastic cover from the tray base, place the air channel adapter (1) into the fraction collector tray base. Make sure the adapter is fully pressed down. This assures that the cold airstream from the thermostat is correctly guided to the tray area of the fraction collector. Place the plug channel (2) on top of the air channel adapter. Both devices must be installed correctly, to assure proper operation of the instrument.

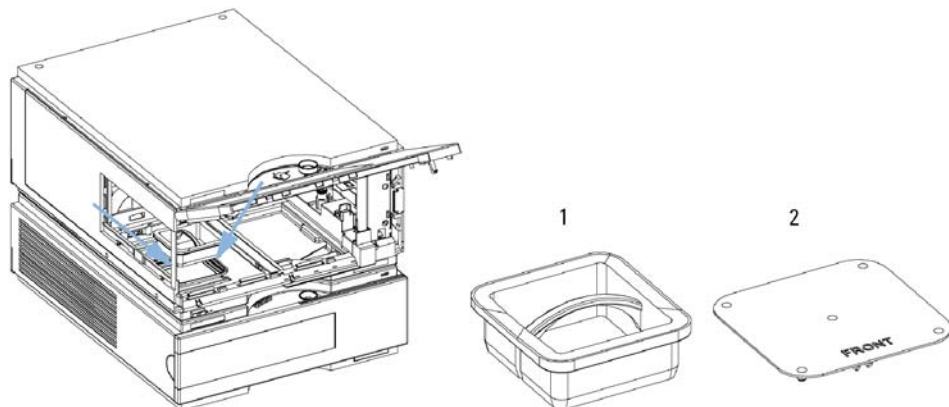


Figure 6 Installation of thermostat and fraction collector

- 11** The fraction collector is delivered with a pre-installed tray compartment divider. This divider should only be used with the thermostatted fraction collector, if a (half or std.) tray is installed into the left and center positions of the instrument. If any tray is installed to the right side of the instrument, remove the tray compartment divider. The tray compartment divider optimizes the cooling performance of the instrument, if only the left and center position of the fraction collector are in use.

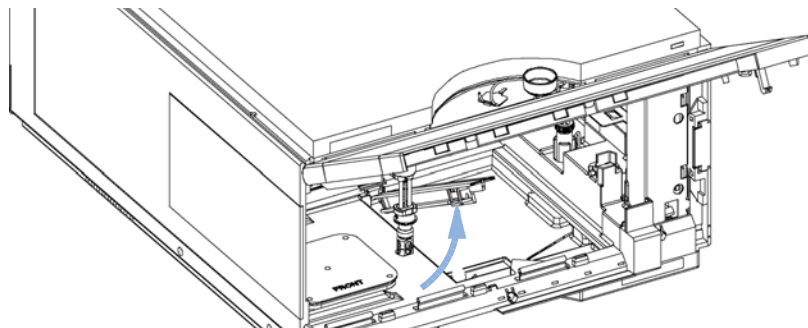


Figure 7 Removing the Tray Compartment Divider

3 Installing the Fraction Collector

Installing a Thermostatted Fraction Collector

- 12** Install the tray you have ordered for your fraction collector. The test tube trays can be adjusted in height, depending on the height of the used test tubes. To adjust the height of the trays, press against the snappers at the guides in the four corners of a tray and move the top plate of the tray up or down (see [Figure 4](#) on page 33).
- 13** Ensure the power switch on the front of the fraction collector is OFF and the power cables are disconnected.
- 14** Connect the cable between the fraction collector and the thermostat, see [Figure 8](#) on page 39.
- 15** Connect the power cables to the power connectors.
- 16** Connect the CAN cable to other Agilent 1260 Infinity modules.
- 17** If an Agilent ChemStation is the controller, connect the LAN connection to the LAN interface.
- 18** Connect the APG remote cable (optional) for non Agilent 1260 Infinity instruments.
- 19** Turn ON power by pushing the button at the lower left hand side of the fraction collector.

- 20** The exhaust fan will turn ON and remove potential solvent vapor from the inside of the instrument. After 2 minutes close the front door. Then the fraction collector will start the hardware initialization process. At the end of this process the status LED should be green.
- 21** The fraction collector is turned ON when the line power switch is pressed and the green indicator lamp is illuminated. The fraction collector is turned OFF when the line power switch is protruding and the green light is OFF.

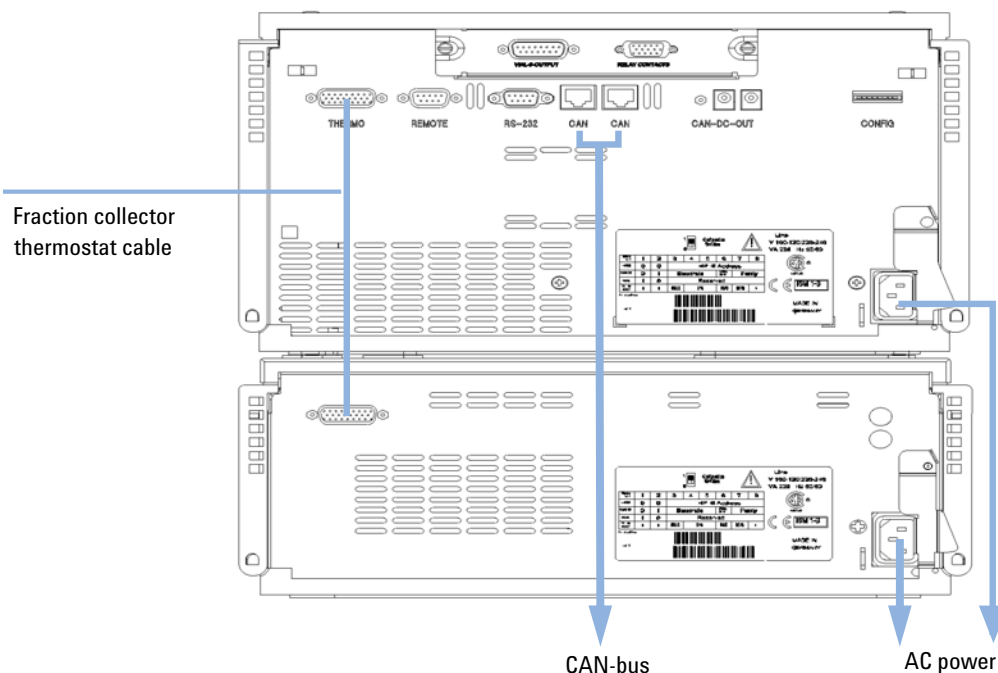


Figure 8 Connections at the Rear of the Thermostatted Fraction Collector

Flow Connections to the Fraction Collector

Parts required	Description Parts from the accessory kit
Preparations	Fraction Collector is installed in the LC system

WARNING

Explosive gas mixtures

There is a risk of explosive gas mixtures in the instrument if flammable solvents are used.

- Cover the plates.
 - Remove the plates from the fraction collector after turning it OFF.
 - Only use solvents with a flash point higher than 200 °C.
-

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.
 - The volume of substances should be reduced to the minimum required for the analysis.
 - Do not operate the instrument in an explosive atmosphere.
-

CAUTION

Liquid spills or fraction losses

Worn or damaged tubings can cause potential spills or lead to fraction losses.

- Explicitly follow the described installation procedures to maximize the lifetime of inlet / waste tubing assembly and the valve to needle tubing, and to avoid potential spills or fraction losses.
- Regularly inspect the tubings and exchange them if they are worn out or show visible signs of damage.

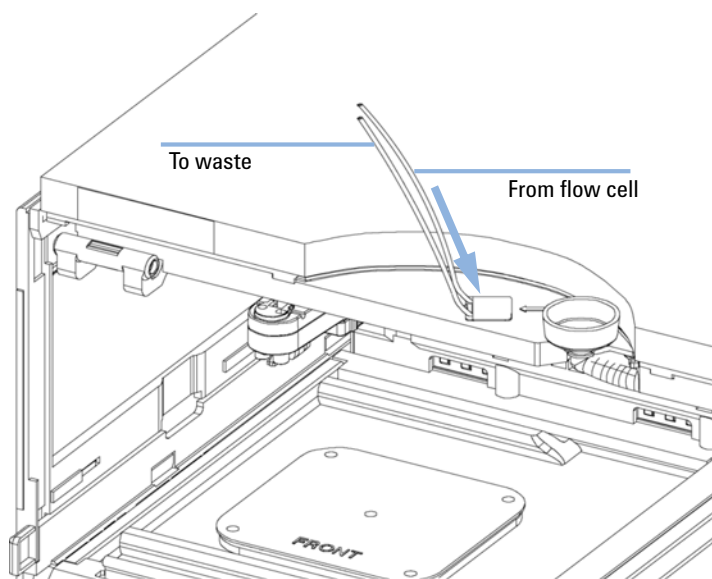


Figure 9 Flow Connections to and from the Fraction Collector

- 1 Route the waste outlet around the instrument as shown below and into a corrugated waste tubing. Fix the waste outlet tubing and the corrugated waste tubing to the instrument with the sticking holders (part of the

3 Installing the Fraction Collector

Flow Connections to the Fraction Collector

Accessory Kit (p/n G1364-68755)) as shown below. Route the corrugated waste tubing into a waste container.

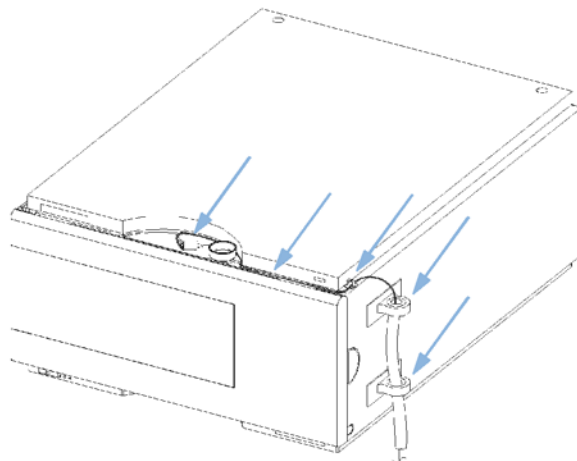


Figure 10 Fixing and Routing the Waste Tubings

Fraction Collector Trays

Installing the Fraction Collector Trays

WARNING

Explosive gas mixtures

There is a risk of explosive gas mixtures in the instrument if flammable solvents are used.

- Cover the plates.
 - Remove the plates from the fraction collector after turning it OFF.
 - Only use solvents with a flash point higher than 200 °C.
-

WARNING

Contamination with adhesives

Closing mats with adhesives can give some contamination in the system. The adhesive is soluble in most of the solvents used in HPLC.

- In general do not use closing mats with adhesive. The fraction collector has no prepunch needle, therefore the adhesive will clog the needle after several injections.
-

- 1** Press the front door latch-holding button located at the front of the right-side cover.
- 2** Lift the front door.
- 3** Adjust the top plate of the test tube tray for the correct tube height if required.
- 4** Load the fraction collector tray with fraction collector well-plates, test tubes or vials as required.
- 5** Slide the fraction collector tray into the fraction collector so that the rear of the tray is seated firmly against the rear of the tray area.

NOTE

Installed trays are automatically detected and identified.

3 Installing the Fraction Collector

Fraction Collector Trays

- 6 Press the front of the fraction collector tray down to secure the tray in the fraction collector.
- 7 Close the front door.

NOTE

If the tray pops out of position the air channel adapter is not correctly inserted.

NOTE

Before starting a run, the instrument has to be correctly configured in the user interface.

Numbering of Vial, Test Tube and Well-plate Positions

With the 4 plates full tray

- Plate in the left front position: P1
- Plate in the left back position: P2
- Plate in the right front position: P3
- Plate in the right back position: P4
- Vessel: A1; A2;... B1; B2;...

With the 2 plates / 10 x 2ml vials or 10 funnels std. trays

- Plate in the front position: P1
- Plate in the back position: P2
- Vessel: A1; A2;... B1; B2;...
- Vials / funnels: 1 - 10

With the 100 vials std. tray

- Vial: 1 - 100

With the half-trays

- Left-hand 40-position tray: 1 - 40
- Center 40-position tray: 101-140
- Right-hand 40-position tray: 201 - 240

or

- Left-hand 15-position tray: 1 - 15
- Center 15-position tray: 101-115
- Right-hand 15-position tray: 201 - 215

With the 40, 60, 125 or 215 position test tube full trays

- Numbering starts in front left corner in columns to the back and then to the right.

Configure Well-plate Types

If the plate you are using is not found on the “[List of Recommended Plates and Closing Mats](#)” on page 171 you may configure a custom plate. Measure the exact dimensions of the plate as marked below and enter the values in the plate configuration table of the ChemStation or the Agilent Instant Pilot G4208A.

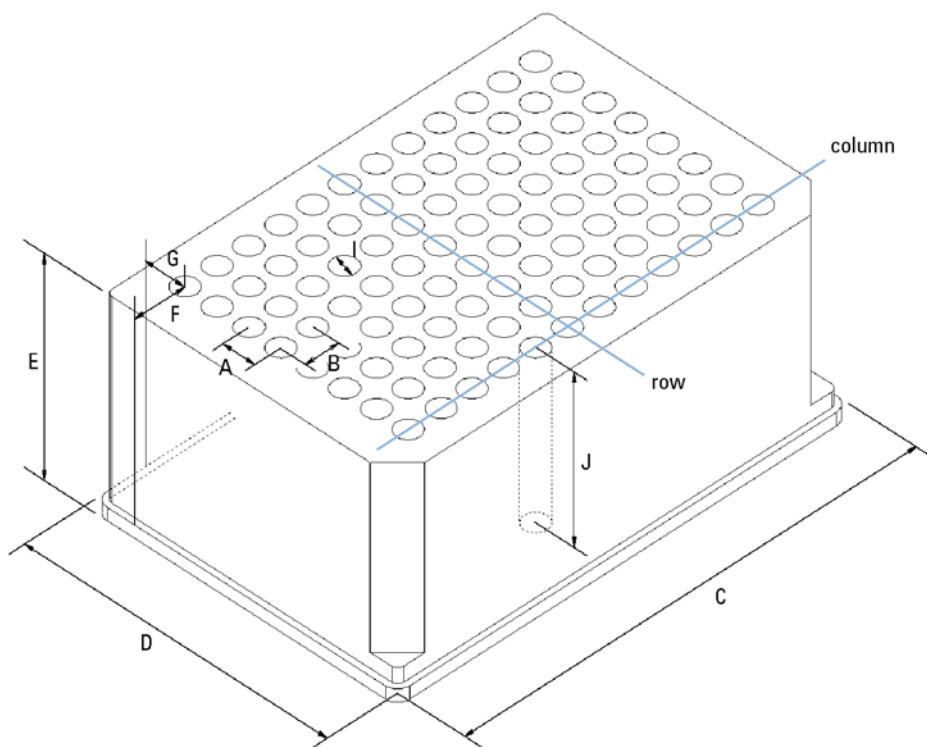


Figure 11 Well-plate Dimensions (straight)

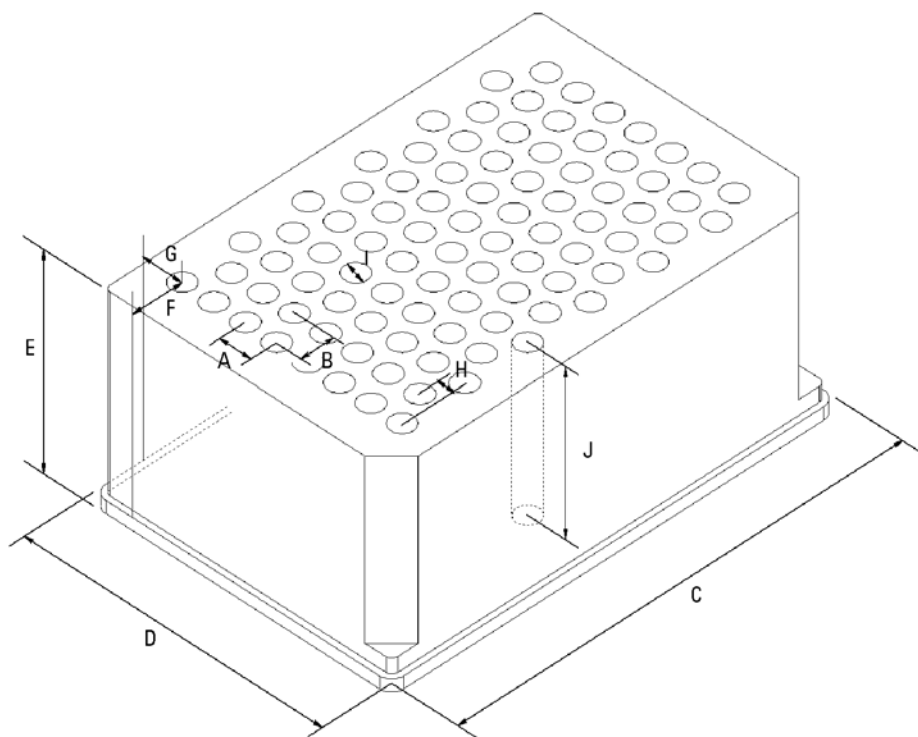


Figure 12 Well-plate Dimensions (staggered)

Table 6 Configuring Well-plate Types

Location	Description	Definition	Limits
	Rows	Number of rows on the plate	up to 16
	Columns	Number of columns on the plate	up to 24
	Volume	Volume (in μl) of a sample vessel	
A	Row distance	Distance (in mm) between the center of two rows	
B	Column distance	Distance (in mm) between the center of two columns	
C	Plate length	X size (in mm) at the bottom of the plate	127.75 $\pm$ 0.25 mm (SBS Standard)

3 Installing the Fraction Collector

Configure Well-plate Types

Table 6 Configuring Well-plate Types

Location	Description	Definition	Limits
D	Plate width	Y size (in mm) at the bottom of the plate	85.50 +/-0.25 mm (SBS Standard)
E	Plate height	Size (in mm) from the bottom to the top of the plate. If you use well plates with inlets, caps and septa, you have to specify the distance from the bottom to the top of the caps.	up to 47 mm
F	Column offset	Distance (in mm) from the left edge (bottom) to the center of the first hole (A1)	
G	Row offset	Distance (in mm) from the back edge (bottom) to the center of the first hole (A1)	
H	Column shift	Offset (in mm) to Y when the rows are not straight but staggered	
J	Well depth	Distance (in mm) from the top of the plate to the bottom of the well. If you use well plates with inlets, caps and septa, you have to specify the distance from the top of the septa to the bottom of the inlets.	up to 45 mm
	Well X size	Size of the well in x direction (Plate length) If you use well plates with inlets, caps and septa, you have to specify the x size of the septa.	min. 3.7 mm min. 3.0 mm with position accuracy alignment (micro scale)
	Well Y size	Size of the well in y direction (Plate width). If you use well plates with inlets, caps and septa, you have to specify the y size of the septa.	down to 3.7 mm min. 3.0 mm with position accuracy alignment (micro scale)>
	Bottom size	For round wells, the relative of the top and bottom of the well	1.0: cylindrical well 0.0: conical well

Table 6 Configuring Well-plate Types

Location	Description	Definition	Limits
	Square	Click in the field to specify whether the well is rectangular or round	Yes: rectangular No: round /oval
	Is well plate	Click in this field to specify if this is a well plate or not. Relevant for continuous flow operation.	Yes: well plate or MALDI Target No: Vial Tray or Eppendorf tray

NOTE

The distances need to be measured with high precision. It is recommended to use a caliper. If possible contact the vendor of non-predefined well plates to obtain these dimensions.

Transporting the Fraction Collector

When moving the fraction collector inside the laboratory, no special precautions are needed. However, if the fraction collector needs to be shipped to another location via carrier, ensure:

- ✓ The transport assembly is in the park position. Use the ChemStation or the Agilent Instant Pilot G4208A for this command.

To move the arm to the park position:

- 1 Switch to the **Diagnosis** view of the ChemStation and select **Fraction Collector > Maintenance Positions** from the **Maintenance** menu.
 - 2 In the upcoming dialog box click **Park Arm**.
- ✓ The vial tray and the sample transport mechanism is secured with the transport protection foam.



4 Using the Fraction Collector

Configuration of the Fraction Collector	52
Delay Volumes and Delay Calibration	58
Detector Signal Delay	59
Performing a Delay Calibration with an UV Detector	62
Setting up a Fraction Collector Method	63
Fraction Preview	67
Starting Your Run with Fraction Collection	69
Viewing Your Results	74
Special Applications	77
Limitations and How to Avoid Problems	80
Check-out Procedures	81
Application Notes	84
Solvent Information	85

This chapter provides information on how to set up the module.



Configuration of the Fraction Collector

Configuration of the Fraction Collector in the ChemStation

In order to setup or change the configuration parameters of your fraction collector select **More Fraction Collector > Configuration** from the **Instrument** menu or right-click on the fraction collector icon in the graphical user interface. In the resulting **Fraction Collector Configuration** dialog box (Figure 13 on page 53) the configuration of the **Trays**, the **Fraction Delay Volumes**, the **Collection Order**, the **Needle Movement** and the **Well-Plates** can be specified.

Trays

In the online mode of the Agilent ChemStation the type of the installed tray is recognized automatically. In the off-line mode an appropriate tray type can be chosen from a dropdown list. For a selection of supported trays see [“Supported Trays for a Fraction Collector”](#) on page 166.

Tube volume [ml]: Volume of the test tube should be specified 10 % less than the maximum fill volume to avoid overflow.

Tube height [mm]: The exact test tube height has to be specified. For the analytical scale fraction collector with 50 mm needle (standard) the maximum height is 48 mm.

NOTE

In the online ChemStation the installed tray is recognized and the default settings for tube volume and tube height are loaded and displayed.

Fraction Collector Configuration : LC113

Trays
 Tray A: 40 Tubes
 Tray B: ---
 Tray C: ---
 Tube Volume [ml]: 45.00
 Tube Height [mm]: 98

Fraction Delay Volumes

	Detector	Volume[µl]
1	DAD1	500

Collection Order
☐ Row by row
☐ Column by column
☒ Shortest path by column
☐ Shortest path by row

Collection Mode
☒ Discrete fractions
☐ Continuous flow

Needle Movement
☐ into location Depth: 5.0 mm
☒ above location Distance: 2.0 mm
☐ contact control Distance: 0.0 mm

Well-Plates
 Plate Type:
☐ Treat as 4x96
 Installed Plates:
☐ Plate 1 (front left) ☐ Plate 3 (front right)
☐ Plate 2 (back left) ☐ Plate 4 (back right)
 Reserved Locations:
 Row(s):
 Column(s):
 Single Location(s):

OK Cancel Help More >>

Figure 13 Fraction Collector Configuration dialog

Use for Sample Recovery: This checkbox only appears, if multiple fraction collectors are configured (see [Figure 24](#) on page 78). Then the last fraction collector can be selected for sample recovery. If four fraction collectors are configured, the fourth one will automatically be used for sample recovery. To change the order of your fraction collectors select **Configure 1260 Access** from the **Instrument** menu. In the **Configuration** dialog box use **Up** or **Down** to change the configuration order and restart your ChemStation.

Fraction Delay Volumes

The **Fraction Delay Volume** specifies the volume between the detector cell and fraction collector diverter valve. This volume has to be determined during the delay calibration. See “[Delay Volumes and Delay Calibration](#)” on page 58 for a detailed description.

Collection Order

The **Collection Order** describes the way of capillary movement during fraction collection. Four different settings are possible:

- **Row by row**
- **Column by column**
- **Shortest path by column**
- **Shortest path by row**

Shortest path by row/column is recommend, if **Continuous flow** is selected as **Collection Mode**.

Collection Mode

The **Collection Mode** defines the triggering of the diverter valve. For **Discrete fractions** the diverter valve switches to waste while the needle moves to the next test tube, vial or well. For **Continuous flow** the diverter valve is not switching during needle movements, except for long movements in the **Row by row** or **Column by column** mode. In addition the needle cannot move **into location** during **Continuous Flow** operation. **Continuous flow** is only available for well plates.

Needle Movement

Into location: In the **Into Location** mode the needle tip moves into the well to the specified **Depth** (in mm). This option is only available for the analytical scale fraction collector in the **Discrete fractions** mode.

Above location: In the **above location** mode the needle tip stays at the specified **Distance** (in mm) above the well during fraction collection.

Contact control: In this mode the needle tip moves down to the well bottom until it reaches the specified **Distance** (in mm) between needle tip and the vial/well bottom. This ensures that the forming droplet is in contact to the vial/well

bottom. During the continuing filling process the needle tip moves upwards while staying in contact with the surface of the collected liquid. This option is recommend for low flow rates and small fraction volumes to avoid air bubbles and accomplish an accurate fraction volume.

Well-Plates

In the **Well-Plates** section the type of well plates used in a well plate tray can be configured. The well plate type can be chosen from the **Plate Type** dropdown list. More detailed information about all pre-configured well plates can be found in the **Instrument** menu. Only one type of well plate can be used on the well plate tray.

The checkbox **Treat as 4 x 96** is available only for 384 well plates and allows to split the 384 well plate virtually into 4 separate 96-well plates. This only changes the collection order. The numbering of the wells remains as indicated on the plate. The four virtual 96 well plates start at locations A1, A13, I1 and I13, respectively.

The filling order of each quarter is as specified in the **Collection Order** section. When the 384 well plate is split into four equal quarters the order of the four plates is the same as displayed in the **Installed Plates** section.

In the **Reserved Locations**, you have the possibility to specify locations that will not be used for Fraction Collection (see [Table 7](#) on page 56).

Table 7 Syntax for the definition of Reserved Locations

Location	Syntax	Description
Rows	A	Row A can't be used
	A, B	Rows A and B can't be used
	A-D	A, B, C and D can't be used
	A-D, F	Rows A, B, C, D and F can't be used
Columns	1	Column 1 can't be used
	1, 2	Columns 1, and 2 can't be used
	1-4	Columns 1, 2, 3 and 4 can't be used
	1-4, 12	Columns 1, 2, 3, 4 and 12 can't be used
Single locations	G12, H12	Locations G12 and H12 can't be used

Configuring Multiple Fraction Collectors

To increase the capacity of the systems up to three fraction collectors can be configured by using the Agilent 1260 Infinity Agilent 12-Position Selector Valve G1160A. In addition a fourth fraction collector can be configured for recovery collection.

- The **Configuration** has to be edited for all fraction collectors. The last fraction collector in the configuration can be selected for sample recovery. To configure the recovery fraction collector please read [“Sample Recovery”](#) on page 78.
- With multiple fraction collectors a fraction start location has to be specified to define the starting position. For further details see [“Assignment of Start Location for Fraction Collection”](#) on page 70. Never use the valve settings to define a start location on systems with multiple fraction collectors.
- We recommend to use inlet tubings of the same length for all fraction collectors. Otherwise a **Delay Calibration** has to be performed for each of those fraction collectors. For the recovery fraction collector the delay volume parameter will be ignored.

NOTE

The **Delay Calibration** can only be performed for fraction collector 1 of your configured system.

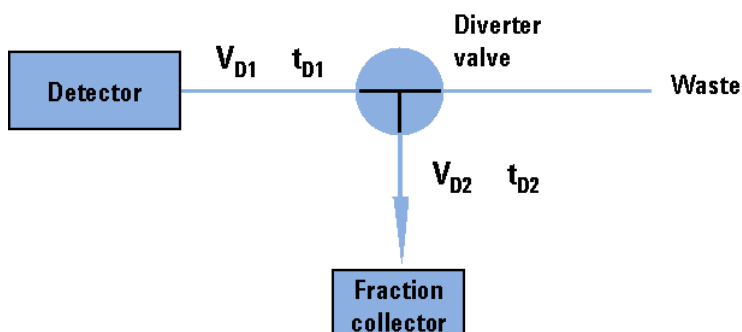
The calculated delay volume is used for all fraction collectors; therefore, the volume (tubing) between the selection valve and each installed fraction collector must be identical.

Otherwise fraction collector 2 and fraction collector 3 have to be temporarily configured as fraction collector 1 during the delay calibration.

-
- The fraction collector inlet tubings have to be connected to port 1-3 of the selection valve in the same order as configured in the ChemStation. The recovery fraction collector is always connected to the waste tubing of the main fraction collectors. The waste tubing from multiple fraction collectors and the inlet tubing to the recovery collector have to be connected through a T-piece.
 - Fraction collector 4 will always be used as a recovery fraction collector.

Delay Volumes and Delay Calibration

Figure on page 58 shows a schematic drawing of the flow path between the detector and the fraction collector with the two delay volumes V_{D1} and V_{D2} . For peak-based fraction collection the system delay times t_{D1} and t_{D2} can be calculated by dividing the delay volumes by the flow rate.



The delay volume V_{D2} is a system parameter, it depends on the installed fraction collector tubing and the fraction collector needle. It is approximately 23 μl . Delay volume V_{D1} , which is specified in the **Fraction Collector Configuration** window, is determined using the **Delay Volume Calibration** feature of the ChemStation software.

When a peak is detected during a purification run (Figure 14 on page 59) the diverter valve is triggered using the following delay time calculations:

- Start of fraction collection: $t = t_0 + t_{D1}$
- End of fraction collection: $t = t_E + t_{D1} + t_{D2}$

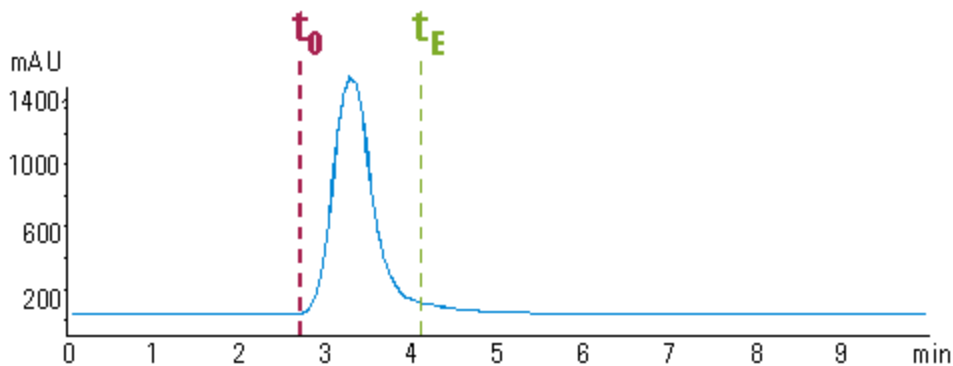


Figure 14 Chromatogram with peak start t_0 and peak end t_E

Detector Signal Delay

Every Agilent 1260 Infinity detector that is used for triggering fractions has an internal signal delay caused by filtering the raw data. The signal delay depends on the **Peakwidth** setting of the detector and is accounted for when the diverter valve is triggered. [Table 8](#) on page 59 to [Table 12](#) on page 61 list the internal signal delay times for different **Peakwidth** settings.

Table 8 Signal Delay Times for the Agilent 1260 Infinity DAD/MWD (G1315D; G4212B/G1365D)

Peakwidth (min)	Response time (sec)	Signal delay (sec)
<0.01	0.1	0.05
>0.01	0.2	0.15
>0.03	0.5	0.5
>0.05	1.0	1.25
>0.10	2.0	2.75
>0.20	4.0	5.9
>0.40	8.0	11.9
>0.85	16.0	23.9

4 Using the Fraction Collector

Delay Volumes and Delay Calibration

Table 9 Signal Delay Times for the Agilent 1260 Infinity DAD/MWD SL (G1315C/G1365C)

Peakwidth (min)	Response time (sec)	Signal delay (sec)
<0.0025	0.02	0.0375
>0.0025	0.05	0.0625
>0.005	0.1	0.144
>0.01	0.2	0.294
>0.03	0.5	0.619
>0.05	1.0	1.27
>0.1	2.0	2.57
>0.2	4.0	5.17
>0.4	8.0	10.4
>0.85	16.0	20.8

Table 10 Signal Delay Times for the Agilent 1260 Infinity VWD G1314B

Peakwidth (min)	Response time (sec)	Signal delay (sec)
<0.005	<0.12	0.07
>0.005	0.12	0.14
>0.01	0.25	0.29
>0.025	0.5	0.58
0.05	1	1.3
0.1	2	2.8
0.2	4	6.0
0.4	8	12.3

Table 11 Signal Delay Times for the Agilent 1260 Infinity VWD SL G1314C

Peakwidth (min)	Response time (sec)	Signal delay (sec)
<0.00125	<0.02	0.0182
>0.00125	0.02	0.0364
>0.0025	0.05	0.0728
>0.005	0.12	0.146
>0.01	0.25	0.328
>0.025	0.5	0.710
>0.05	1	1.49
>0.1	2	3.08
>0.2	4	6.26
>0.4	8	12.6

Table 12 Signal Delay Times for the Agilent 1260 Infinity FLD G1321A

Peakwidth (min)	Response time (sec)	Signal delay (sec)
<0.005	<0.12	0.018
>0.005	0.12	0.073
>0.01	0.25	0.18
>0.03	0.5	0.44
>0.05	1	0.96
>0.1	2	2.0
>0.2	4	4.2
>0.4	8	8.6

4 Using the Fraction Collector

Delay Volumes and Delay Calibration

CAUTION

Loss of data

If the internal signal delay is longer than the delay time t_{D1} some of the peak will be lost.

→ The maximum allowed signal delay time can be calculated using the following equation:

$$\text{Signal delay time}_{(max)} = \frac{V_{D1}}{\dot{v}} \quad \dot{v} = \text{Flow rate}$$

→ After calculating the maximum signal delay time a **Peakwidth** setting can be selected that gives a signal delay time, which is shorter than the calculated maximum signal delay time. This **Peakwidth** setting should then be used for the LC purification run.

NOTE

We recommend to set the **Peakwidth** always to > 0.01 for the DAD and MWD or to > 0.005 for the VWD.

If the **Peakwidth** setting cannot be reduced and the signal delay time is longer than t_{D1} it is also possible to increase V_{D1} by adding additional tubing. However this higher delay volume will increase the peak dispersion between detector and fraction collector.

The stop-time of the run in the ChemStation must be set to at least:

Total duration of time table (time of last entry OFF) + fraction collector delay time ($V_{D1}/\dot{v}$) + 0.1 min for time-based fraction collection.

End of last peak (t_E) + fraction collector delay time ($V_{D1}/\dot{v}$) + 0.1 min for peak-based fraction collection

Performing a Delay Calibration with an UV Detector

- 1 Place a vial containing the Delay sensor calibrant (p/n G1946-85020) in position 1 of the autosampler.
- 2 Perform Delay Calibration by using Lab Advisor B.02.01 or higher. Follow instruction given in Lab Advisor.

Setting up a Fraction Collector Method

In order to setup or change the method parameters of your fraction collector select **Setup Fraction Collector** from the **Instrument** menu or right-click on the fraction collector icon in the graphical user interface. This will open the **Setup Fraction Collector** dialog box as displayed in [Figure 15](#) on page 63. In the **Setup Fraction Collector** dialog box general method settings are specified.

Set up Fraction Collector : LC113

Fraction Trigger Mode

☐ Off ☐ Use Timetable ☒ Peak-based Max. Peak Duration min

Peak Detectors

	Detector	Working Mode	Up Slope [Unit/s]	Down Slope [Unit/s]	Threshold [Unit]	Upper Threshold [Unit]	Unit
1	DAD1	Threshold/Slope	5.00	10.00	10.000	3000.000	mAU

Fraction is collected when a peak is detected by:

☐ all selected peak detectors ☒ at least one selected peak detector ☐ Use MSD for Mass-based Fraction Collection

Timetable

Time	Trigger Mode	Max. Peak Duration [min]	# of Fractions	Timeslices [min]
------	--------------	--------------------------	----------------	------------------

Time

Stoptime: as Pump min
Posttime: Off min

Rinse Fraction Collection Needle

☐ At the start of collection
☐ Between fraction collection

Auxiliary

Max. fill volume per location as configured ml

☐ Store temperature signal
☐ Store delay sensor signal
☐ Store UJB signal

Figure 15 Set up Fraction Collector dialog box

Fraction Trigger Mode

Use Timetable: Enables the **Timetable**.

Peak-based: If **Peak-based** is selected, the collection of a fraction is triggered by the signal of the detector, e.g. the Agilent 1260 Infinity Diode Array Detector or Variable Wavelength Detector. The detailed trigger conditions are specified in the **Peak Detectors** table. In the peak-based trigger mode all entries in the timetable are ignored.

Max. Peak Duration: Defines a maximum collection time in case that the signal does not reach the condition to cut the fraction as exhibited in [Figure 16](#) on page 64. This could be caused by tailing peaks or if the baseline is drifting during gradient runs. The default value is set to 0.5 minutes. If broad peaks are expected, this value should be increased without exceeding the run time.

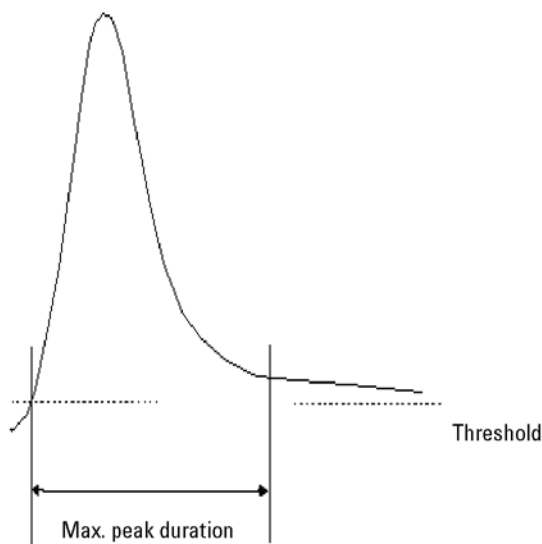


Figure 16 Maximum Fraction Duration

Peak Detectors

In the **Peak Detectors** section a list of all peak detectors that are connected to the system is displayed. Agilent 1260 Infinity diode array detectors, multiwavelength detectors, variable wavelength detectors and fluorescence detectors are recognized automatically. Other detectors, e.g. Agilent 6000 mass-selective detectors or HP1050 detectors, are connected through the Universal Interface Box (UIB).

The peak detector table contains seven columns:

Working Mode

For each peak detector **Threshold only**, **Threshold/Slope** or **Slope only** are possible.

In the **Threshold only** mode the settings for **Up Slope**, **Down Slope** and **Upper Threshold** in the subsequent columns are ignored. Fraction collection is triggered whenever the detector signal exceeds the specified threshold value. When the signal drops below the threshold value fraction collection is stopped.

In the **Slope only** mode fraction collection is triggered on the slope of the detector signal. Adequate values for **Up Slope** and **Down Slope** can be specified in the corresponding fields.

In the **Threshold/Slope** mode fraction collection is triggered on the corresponding values for threshold and slope. The fraction collection is started if the detector signal exceeds both the threshold and the **Up Slope** value. The fraction collection is stopped if the detector signal drops either below the threshold or the **Down Slope** value.

To specify the trigger values **Up Slope**, **Down Slope**, **Threshold** and **Upper Threshold** we recommend to use the **Fraction Preview** tool as described in “[Fraction Preview](#)” on page 67.

Upper Threshold

At high absorbance values the light intensity on the detector is extremely low and consequently detector noise will be superimposed on the detector signal. In this case the detector noise might trigger fraction collection. To avoid false fraction collection triggering, we recommend setting an **Upper Threshold** well below the limit where this false triggering effect might occur. As soon as the detector signal exceeds the **Upper Threshold**, settings for **Up Slope** or **Down Slope** will be ignored until the signal drops again below the **Upper Threshold**.

4 Using the Fraction Collector

Setting up a Fraction Collector Method

When using more than one peak detector fraction collection can be triggered either when **all selected peak detectors** detect a peak or when **at least one selected peak detector** detects a peak basing on the settings in the peak detectors table above.

If an MSD is used for mass-based fraction collection, **Use MSD for mass-based Fraction Collection** must be checked.

Timetable

The **Timetable** can be used to program changes in the Fraction Trigger Mode during the analysis by entering a Time and specifying the trigger settings.

Trigger Mode Off, Peak Based and Time Based can be selected. If the Off is selected, no fractions are collected. The last entry in the timetable has to be the command Off.

Whenever the **Peak Based** mode is specified fractions will be collected based on the peak detection parameters given in the Peak Detector table. Additionally a **Maximum Peak Duration** in minutes has to be specified. This parameter is mandatory if you use Peak Controlled fraction collection, but is disabled for Time Based fraction collection.

When the Time Based mode is chosen two different options are available:

- The **# of Fractions** can be edited to collect a fixed number of equal fractions in a give time interval. This time interval is defined by the time value in the current and following timetable line.
- **Timeslices [min]** can be edited to collect fractions with a defined collection time. With this option the collection time of the last fraction can be shorter. This depends on the overall runtime.

For editing the Timetable the functions **Insert**, **Append**, **Cut**, **Copy** and **Paste** are offered.

To access the additional sections in the **Setup Fraction Collector** dialog box click **More**.

Time

In the time section of the dialog box the **Stoptime** and the **Posttime** for the fraction collector can be specified. By default the Stoptime is set to as pump and the posttime is switched OFF.

Rinse Fraction Collection Needle (Analytical Scale only)

If **Discrete fractions** is selected as **Collection Mode** (see also “[Collection Mode](#)” on page 54), you can setup a needle rinse step before the fraction collection and/or between fractions. Then the needle will move to the funnel on the internal tray and the diverter valve will switch to flush the needle in order to avoid carry over from the previous fraction. The instrument will determine, if it is possible to rinse the needle before the next fraction is expected.

If you have recovery positions in your fraction collector or if you are using one fraction collector for sample recovery in a multiple fraction collector configuration, the function **rinse between fractions** is ignored.

Auxiliary

In the Auxiliary section the **Maximum fill volume** per location can be specified. If as configured is selected, the pre-configured volume (see **Instrument > Pre-configured Wellplate Types**) is used. This ensures that the location (well, vial or tube) cannot be overfilled during fraction collection. This volume can be further reduced by defining a customized volume.

Additional check boxes in this section provide the opportunity to **Store temperature signal** and to **Store UIB signal**.

Fraction Preview

To determine the appropriate fraction collection parameters the Agilent ChemStation provides a valuable tool that becomes accessible by pushing the button labelled Fraction Preview Tool ([Figure 17](#) on page 68) in the Peak Detectors section.

4 Using the Fraction Collector

Setting up a Fraction Collector Method

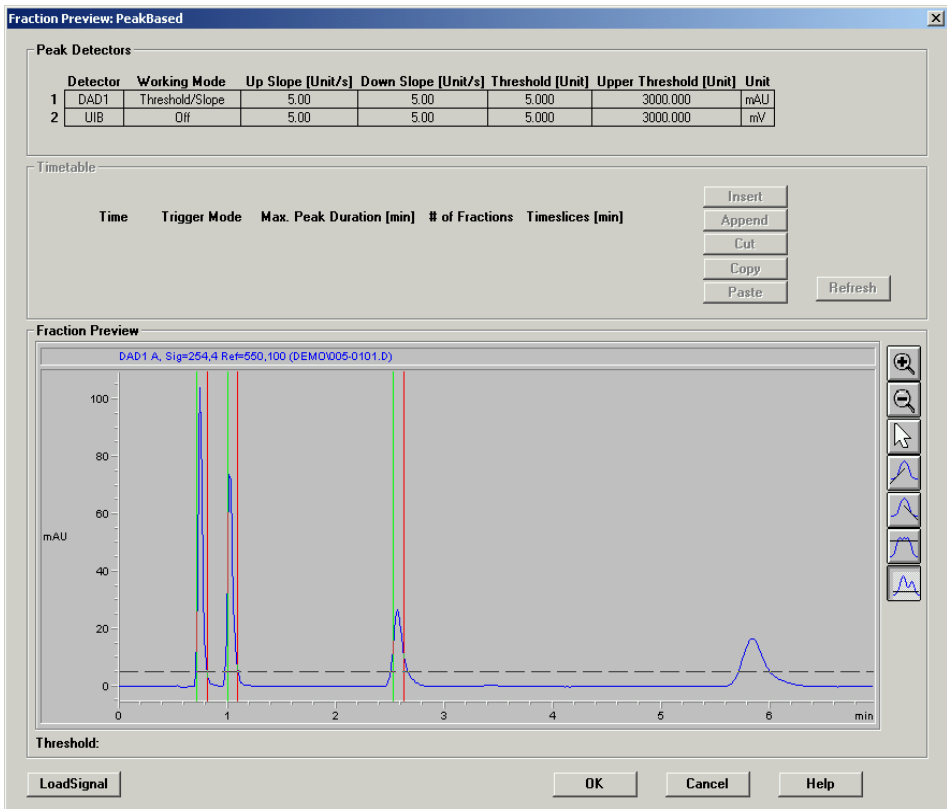


Figure 17 Fraction Preview dialog box

The **Fraction Preview** screen allows to test the fraction collection parameters against an example chromatogram. It can also be used to optimize the fraction collection parameters interactively. With the help of this tool values for up and down slope as well as for upper and lower threshold can easily be graphically specified. To load a chromatogram e.g. a pilot run click **Load Signal**. Parameters can now be changed either manually in the detector table and **Timetable** or graphically in the **Fraction Preview** screen. By clicking the desired buttons on the right hand side of the **Fraction Preview** screen the chromatogram can be zoomed, the values for up and down slope can be specified and the upper and lower threshold level can be set-up. The graphically specified values are automatically transferred to the **Peak Detector** table.

Starting Your Run with Fraction Collection

Resetting the Current Fill Levels

To protect the collected fractions against contaminations and to avoid overfilling of single positions, the fraction collector stores current fill levels. It will block used positions, if no fraction start location has been specified.

If a new tray of the same type is inserted in the fraction collector, all fill levels can be reset (not just the well plate, vial or test tube). After re-installation of the tray a pop-up window (see [Figure 18](#) on page 69) will be displayed. Click **Yes** to reset all fill levels from the previous run.



Figure 18 Reset fill volumes

CAUTION

Contaminated fractions, unwanted overfill events or split fractions

The window ([Figure 18](#) on page 69) to reset current fill levels is not displayed, if the well plate, the vial or the test tube is removed without removing the complete tray. Consequently contaminated fractions, unwanted overfill events or split fractions might be observed during the following LC run.

→ Remove the complete tray to correctly reset fill levels.

Assignment of Start Location for Fraction Collection

The start location for fraction collection can either be assigned in the **Sample Info** dialog box ([Figure 19](#) on page 71) in the **Run Control** menu, in the **Sequence Parameters** dialog box ([Figure 20](#) on page 72) or in the **Sequence Table...** in the **Sequence** menu. All fraction start locations are entered in the following formats.

For vials and test tubes:

fraction collector # - vessel number, e.g. *1-Vial 1*

For well plates:

fraction collector # - plate number - row - column, e.g. *1-P1-A-01*

with

fraction collector # = 1,2; plate number = P1-P4; row is A to H and column = 1-12 for 96-well plates.

NOTE

Note the orientation of the well plate. The starting position A1 is always in the back left corner of the plate.

Sample Info: Instrument 1

Operator Name: Bio Chemist

Data File

☐ Prefix/Counter ☒ Manual

Filename: DEFAULT.D

Subdirectory: PROTEOME

Path: C:\HPCHEM\1\DATA\

Sample Parameters

Location: Vial 1 (blank run if no entry)

Fraction Start Location: 1-P1-A-01

Sample Name: Sample_01 Plate ID:

Sample Amount: 0 Multiplier: 1

ISTD Amount: 0 Dilution: 1

Comment:

Run Method OK Cancel Help

Figure 19 Sample Info dialog box

All settings made in the **Sequence Table** override the settings made in the **Sequence Parameters** screen. In addition to the exact position for fraction collection start in the **Sequence Table** it is also possible to specify **Next Plate** and **Next Location**. In the former case fraction collection is started at the next free plate and in the latter case fraction collection is started at the next free location.

If a fraction start location is specified, the collector will start at this location, even if this vessel has been used during a previous run.

If no fraction start location has been specified, the fraction collection starts at the first empty vessel or location.

4 Using the Fraction Collector

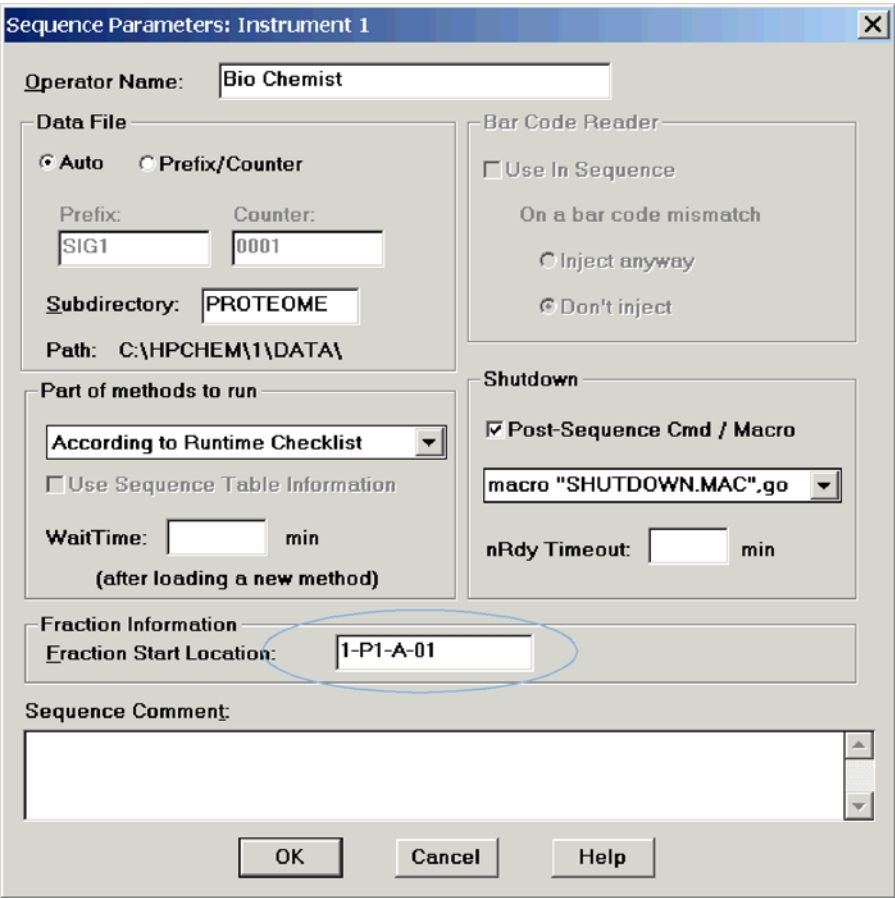
Starting Your Run with Fraction Collection

CAUTION

Loss of data, loss of fractions

With multiple fraction collectors a fraction start location must be specified. Otherwise the fraction collection of the next run will start at the fraction collector, where the last run stopped.

- Always specify a fraction start location, e.g. 1-P1-A-01 or 2-Vial 10.
- Never use the valve settings to define a start location.



The image shows a screenshot of the 'Sequence Parameters: Instrument 1' dialog box. The 'Operator Name' field is filled with 'Bio Chemist'. Under 'Data File', 'Auto' is selected, and the 'Prefix' is 'SIG1' and 'Counter' is '0001'. The 'Subdirectory' is 'PROTEOME' and the 'Path' is 'C:\HPCHEM\1\DATA\'. Under 'Part of methods to run', 'According to Runtime Checklist' is selected in the dropdown, and 'Use Sequence Table Information' is unchecked. The 'WaitTime' is set to 0 minutes. Under 'Fraction Information', the 'Fraction Start Location' is '1-P1-A-01', which is circled in blue. The 'Sequence Comment' field is empty. At the bottom are 'OK', 'Cancel', and 'Help' buttons.

Sequence Parameters: Instrument 1

Operator Name: Bio Chemist

Data File

☒ Auto ☐ Prefix/Counter

Prefix: SIG1 Counter: 0001

Subdirectory: PROTEOME

Path: C:\HPCHEM\1\DATA\

Bar Code Reader

☐ Use In Sequence

On a bar code mismatch

☐ Inject anyway

☒ Don't inject

Part of methods to run

According to Runtime Checklist

☐ Use Sequence Table Information

WaitTime: 0 min

(after loading a new method)

Shutdown

☒ Post-Sequence Cmd / Macro

macro "SHUTDOWN.MAC",go

nRdy Timeout: 0 min

Fraction Information

Fraction Start Location: 1-P1-A-01

Sequence Comment:

OK Cancel Help

Figure 20 Start location in **Sequence Parameters** dialog box

To start a single run

- click **Start** in the graphical user interface of the ChemStation.
- select **Run Method** from the **RunControl** menu.
- press **F5**.

To start the sequence

- click **Start** in the graphical user interface of the ChemStation.
- select **Run Sequence** from the **RunControl** menu.
- press **F6**.

Online Tick Marks

To display tick marks for fraction start and stop events in you Online Plot click **Change** in your **Online Plot** window. Then check **Show fraction collection ticks** in the **Edit Signal Plot** window (Figure 21 on page 73).

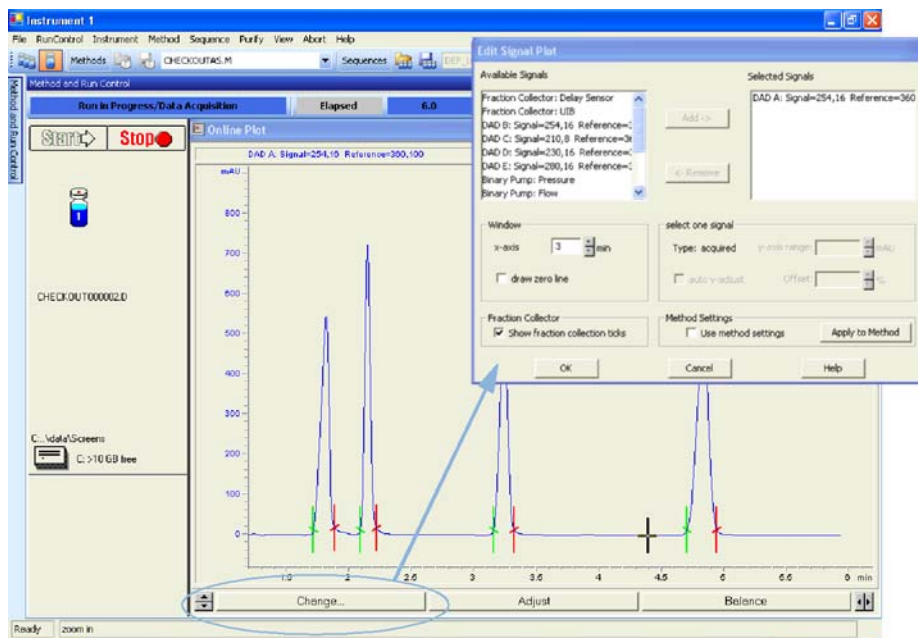


Figure 21 Online tick marks

Viewing Your Results

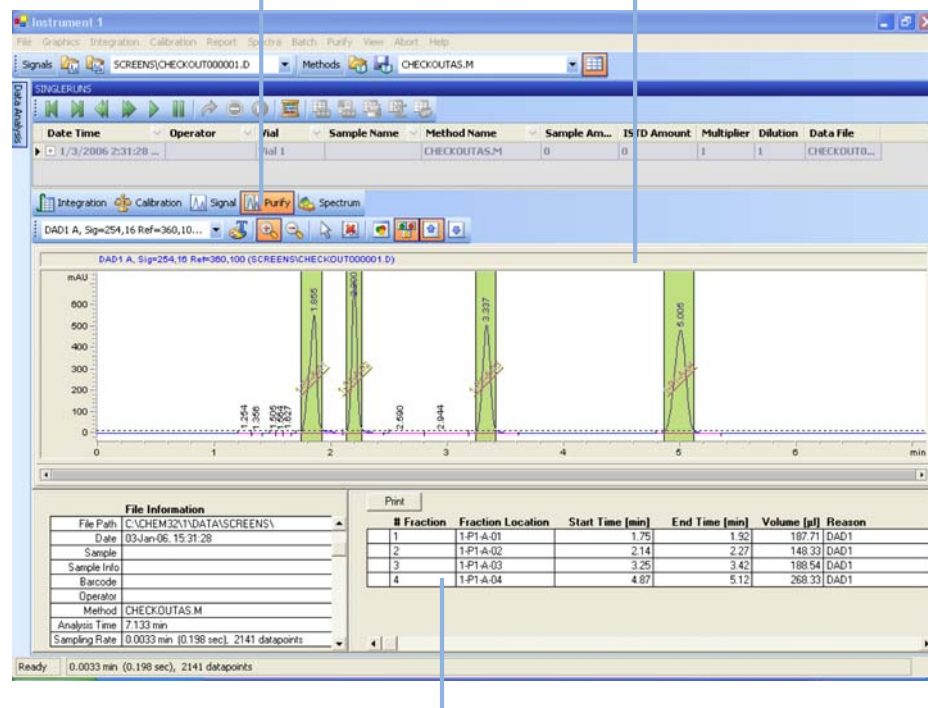
Data Analysis

In order to display the tick marks for the collected fractions on the screen, click **Signal options** from the **Graphics** menu. Then choose **Separated** in the **Layout dropdown list**.

To review your chromatograms, file information and a fraction list, select the **Data Analysis** view and click **Fraction Task** as displayed below.

Chromatogram with Fraction Tick Marks

Purification Task



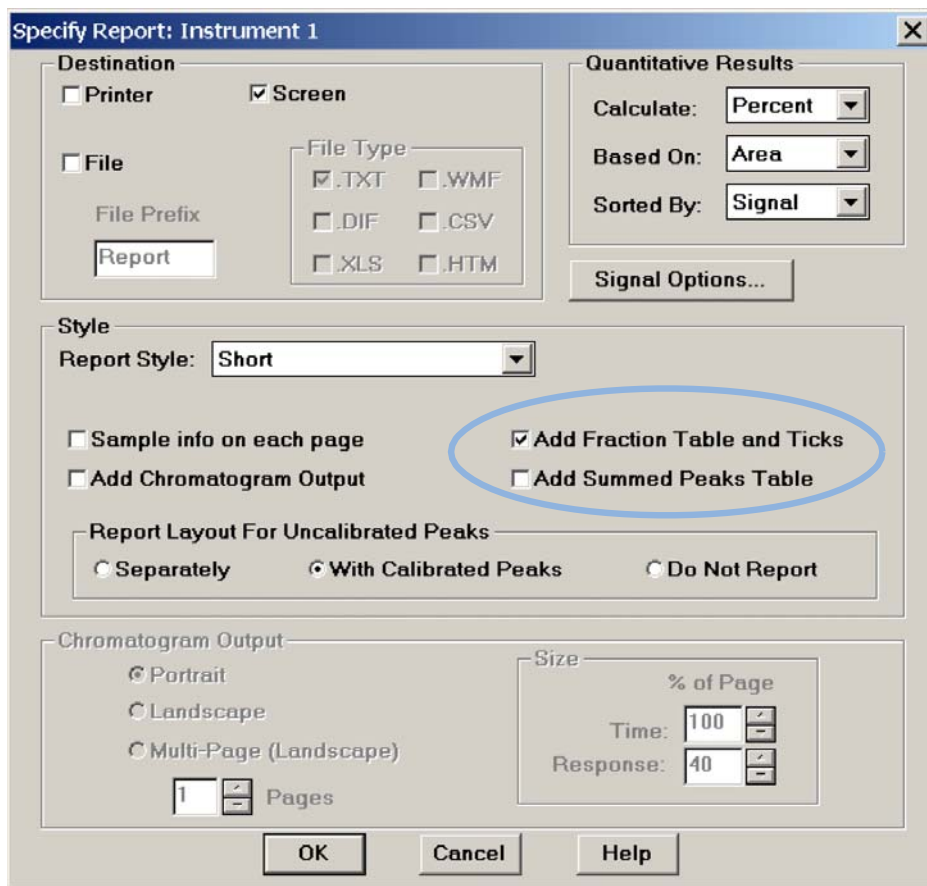
Fraction table

Figure 22 Fraction Task button

In order to display the tick marks for the collected fractions on the screen, click **Signal options** from the **Graphics** menu. Then choose **Separated** in the Layout dropdown list.

Report

In order to create reports with a fraction table and tick marks the item **Add Fraction Table and Ticks** in the **Specify Report** box has to be checked.



The image shows a software dialog box titled "Specify Report: Instrument 1". It contains several sections for configuring report output:

- Destination:** Includes checkboxes for "Printer" and "Screen" (checked). There is also a "File" checkbox and a "File Prefix" text field containing "Report".
- File Type:** A group box containing checkboxes for ".TXT" (checked), ".WMF", ".DIF", ".CSV", ".XLS", and ".HTM".
- Quantitative Results:** Includes dropdown menus for "Calculate:" (set to "Percent"), "Based On:" (set to "Area"), and "Sorted By:" (set to "Signal"). A "Signal Options..." button is located below.
- Style:** Includes a "Report Style:" dropdown set to "Short".
- Content Options:** Includes checkboxes for "Sample info on each page", "Add Chromatogram Output", and "Add Fraction Table and Ticks" (which is circled in blue). There is also an unchecked checkbox for "Add Summed Peaks Table".
- Report Layout For Uncalibrated Peaks:** Includes radio buttons for "Separately", "With Calibrated Peaks" (selected), and "Do Not Report".
- Chromatogram Output:** Includes radio buttons for "Portrait" (selected), "Landscape", and "Multi-Page (Landscape)". Below these is a "Pages" spinner set to "1".
- Size:** A sub-section with the label "% of Page" containing spinners for "Time:" (set to 100) and "Response:" (set to 40).

At the bottom of the dialog are "OK", "Cancel", and "Help" buttons.

Special Applications

Pooling

We define pooling as collecting fractions from multiple injections into the same fraction locations. In the Agilent ChemStation, there are two ways to initiate pooling:

- If you want to collect multiple injections from the one sample vial or sample well into the same fraction location, specify multiple injections in the **Sequence Table**.
- In case you want to collect multiple injections from the several sample vials or sample wells into the same fraction location, specify the same **Fract. Start** for multiple samples in the **Sequence Table**. An example is displayed in [Figure 23](#) on page 77. In this sequence the sample 1 will be injected from plate 1 position A1 of the Agilent 1260 Infinity High Performance Autosampler and fraction collection will start at plate 1 position A1 of the fraction collector.

For the next sample 2, which is injected from a different location in the high performance autosampler, the fraction collection will start again at plate 1 position A1 of the fraction collector.

Line	Location	Sample Name	Dilution	Datafile	Inj Volume	Fract. Start	LimsID	Target Masses
1	P1-A-01	Sample 1				1-P1-A-01		
2	P1-A-02	Sample 2				1-P1-A-01		
3	P1-A-03	Sample 3				1-P1-A-01		
4								

Figure 23 Pooling sequence

CAUTION

Data loss due to pump shut-down

If you specify multiple injections for pooling, the overflow protection is not activated. If a fraction collection location is overflowed, an error message occurs and the pump is shut OFF.

- Make sure that all fraction collection locations are large enough to completely collect all pooled fractions.

Sample Recovery

The Agilent 1260 Infinity Fraction Collector offers different possibilities for sample recovery:

- The preferred recovery strategy is to install multiple fraction collectors in your LC systems and use the last of those fraction collectors for recovery. This recovery fraction collector can be selected in the Configuration dialog box as displayed in [Figure 24](#) on page 78. The fourth fraction collector in your systems will always be used for sample recovery. Also read [“Configuring Multiple Fraction Collectors”](#) on page 56.

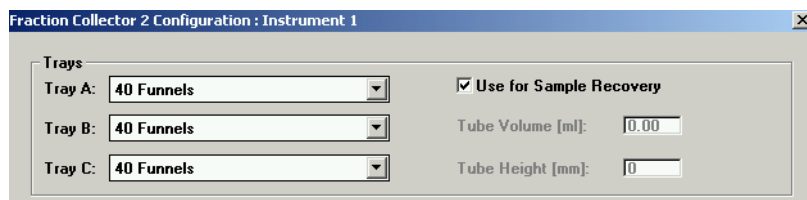


Figure 24 Sample Recovery Fraction Collector

- For the following tray configurations fixed recovery location will be assigned from the ChemStation. In order to disable the recovery the tray configuration has to be changed.
 - a** Standard tray for two well plates + 10 x 2 ml vials (p/n G2258-60011) and Halftray for 40 x 2 mL vials (p/n G1313-44512).
In this configuration the forty 2 ml vials on the half will automatically be used for recovery.
 - b** Standard tray for two well plates + 10 x 2 ml vials (p/n G2258-60011) and Halftray for 15 x 6 mL vials (p/n G1313-44513).
In this configuration the fifteen 6 mL vials on the half will automatically be used for recovery.
 - c** Standard tray for two well plates + 10 x 2 ml vials (p/n G2258-60011) and Half Tray for 40 Funnels (p/n G1364-84532).
In this configuration the funnels on the half will automatically be used for recovery.

- The Std. tray for 2 well plates + 10 collecting funnels (p/n G1364-84522) allows two modes of operation.
 - a** If well-plates are configured in the fraction collector configuration the (checkbox **Plate 1** or **Plate 2** are selected), the 10 funnel locations will automatically be used for recovery.
 - b** If the checkbox **Plate 1** and **Plate 2** are not selected, the 10 funnel position can be used for regular fraction collection.

CAUTION

Fraction contamination

With each start of a new sequence the recovery will start at the same position. This can lead to fraction contamination, if the vessels containing the recovery fractions are not exchanged.

→ Exchange the vessels containing the recovery fractions before starting a new sequence.

NOTE

The number of recovery locations automatically defines the maximum number of injections. When using the standard tray for two well plates and 10 funnels, only ten injections per sequence are executed.

Limitations and How to Avoid Problems

Rinse Fraction Collection Needle

If *Rinse Fraction Collection Needle* is set to *Between fraction collection*, at least 0.3 min are required to perform this task.

When doing time-based fraction collection rinsing the needle is only possible between two time table entries, which must have a gap of at least 0.3 min. For peak-based fraction collection a time gap of also at least 0.3 min is required. If a new peak is detected during the rinse process, it is aborted and the needle moves back to the next free fraction position. Depending on flow rate and delay volume VD1 the beginning of this peak may be lost.

If you have recovery positions in your fraction collector or if you are using one fraction collector for sample recovery in a multiple fraction collector configuration, the rinse function **between fraction collection** is ignored.

Needle Movement

The option **into location** under **Needle Movement** in the fraction collector configuration must only be used for capped 2 or 5 mL vials or well-plates. Using other or open vials with this command can lead to a **Movement failed** error.

Replacing fraction containers

When replacing filled tubes, vials or well-plates from the fraction collector make sure to remove and re-insert the complete tray. Otherwise the fraction collector will not recognize that the fraction containers were emptied.

Pooling

When pooling fractions, overflow protection no longer exists. It is the user's responsibility to make sure that all fraction collection locations are large enough to completely collect all pooled fractions. If a fraction collection location is overfilled, an error message occurs and the pump is shut OFF.

Check-out Procedures

The purpose of the check-out injection is the verification of the correct installation and configuration of the Agilent 1260 Infinity Purification System. In addition the checkout procedure can be used for troubleshooting, if the Agilent 1260 Infinity Fraction Collector doesn't operate as expected.

The checkout procedure depends on the flow range and on the peak trigger (UV-based or mass-based).

In the following sections the check-out procedures for the UV based trigger are described. For the mass-based system refer to the Agilent 1260 Infinity Purification System User's Guide (p/n G2262-90011).

Check-out Procedure for UV-triggered Fraction Collection AS

Prerequisites

- ✓ The Agilent 1260 Infinity LC has been flushed thoroughly with the liquid phase as described in [Table 13](#) on page 82 and the lamp(s) of the UV detector have been switched on for at least 15 minutes to provide a stable baseline.
- ✓ The Agilent 1260 Infinity Fraction Collector has been installed and configured. The delay volume has been calibrated as described in [“Performing a Delay Calibration with an UV Detector”](#) on page 62
- ✓ The 40-vial tray (Halftray for 40 x 2 mL vials (p/n G1313-44512)) with at least four 2 ml vials is installed.
- ✓ All wells are empty and the fill levels, which are stored on the module, have been reset. To verify this remove and insert the 40-vial tray. Then click **Yes** in the displayed dialog box.

Table 13 Method Parameters for UV-triggered Fraction Collection Check-out AS

Agilent 1260 Infinity modules	• Standard or micro degasser • Isocratic, quaternary or binary pump • Autosampler/ well plate sampler • Column compartment • Diode-array detector, multi-wavelength detector or variable wavelength detector with analytical flow cell • Fraction collector analytical scale
Column	Zorbax Eclipse XDB C18, 150 mm x 4.6 mm, 5 µm (p/n 993967-906) or equivalent
Sample	Agilent isocratic checkout sample (p/n 01080-68704)
Flow	0.8 mL/min
Stoptime	6 minutes
Solvent A	10 % (HPLC grade water)
Solvent B	90 % (HPLC grade acetonitrile)
Wavelength VWD	254 nm
Wavelength DAD/MWD	Signal A: 254/16 nm, Reference: 360/80 nm
Peakwidth (Responsetime)	> 0.1 min (2 s)
Injection volume	1 µL
Column temperature	40.0 °C
Fraction trigger mode	Peak-based
Max. peak duration	0.5 min
DAD/MWD working mode	Threshold/slope
Upslope	5 mAU/s
Downslope	10 mAU/s
Threshold/Upper threshold	10 mAU/3000 mAU

Expected Result

Run a single injection under the conditions as described in Table 13 on page 82. You should observe a chromatogram with four peaks as displayed in Figure 25 on page 83. Four fractions should have been collected in the fraction collector vials.

NOTE

The retention times of the peaks could be different depending on the system configuration (e.g. flow cell type, capillary volumes, type and age of column).

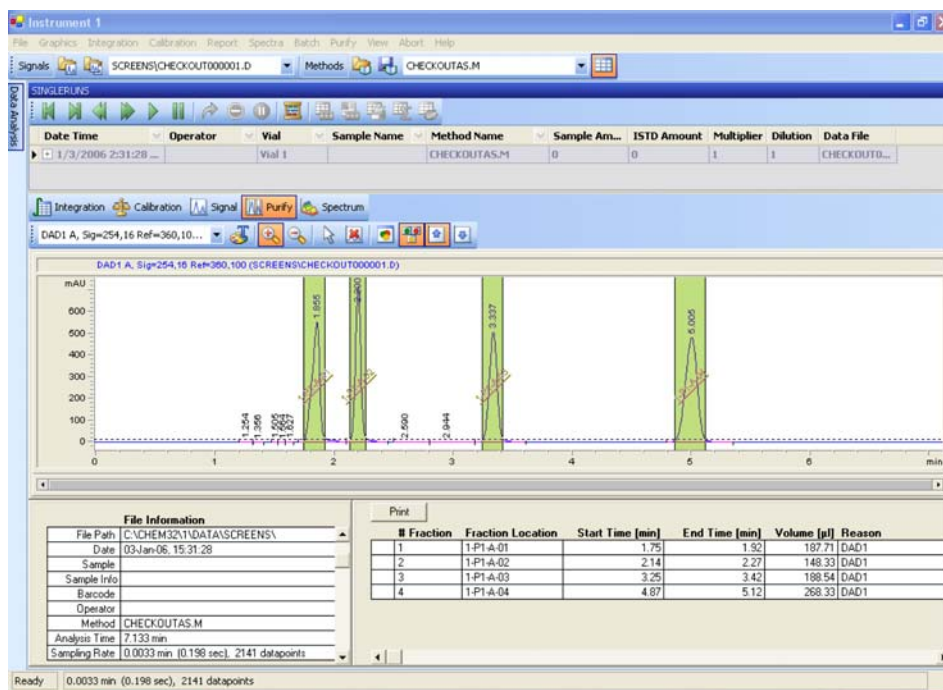


Figure 25 Chromatogram of the check-out injection with UV-based trigger

Application Notes

More information about the Agilent 1260 Infinity fraction collectors and purification systems are available from a of application notes. Printed versions can be ordered from Agilent or pdf-files can be downloaded from the Agilent Website

<http://www.chem.agilent.com>

Table 14 Selected Agilent Technologies Application Notes

Title	Publication Number
An optimized system for analytical and preparative work	5988-9649EN
Recovery collection with the Agilent LC purification system	5988-9650EN
Solutions for preparative HPLC-Application Compendium	5988-9646EN
Innovative fraction collection with the Agilent LC purification platform	5988-9250EN
Automated fraction re-analysis - does it really make sense?	5988-8653EN
Injection of high-concentration samples with the Agilent LC purification system	5988-8654EN
Sophisticated peak-based fraction collection - working with up and down slope	5988-7895EN
Strategies for purification of compounds from non-baseline separated peaks	5988-7460EN
Method scale-up from analytical to preparative scale with the Agilent LC purification system PS	5988-6979EN
Peak-based fraction collection with the Agilent LC purification system AS - Influence of delay volume on recovery	5988-5747EN

Solvent Information

Observe the following recommendations on the use of solvents.

- Follow recommendations for avoiding the growth of algae, see pump manuals.
- Small particles can permanently block capillaries and valves. Therefore, always filter solvents through 0.4 µm filters.
- Avoid or minimize the use of solvents that may corrode parts in the flow path. Consider specifications for the pH range given for different materials like flow cells, valve materials etc. and recommendations in subsequent sections.

Solvent information for parts of the 1260 Infinity Bio-inert LC system

For the Agilent 1260 Infinity Bio-inert LC system, Agilent Technologies uses highest quality materials (see “[Bio-inert Materials](#)” on page 10) in the flow path (also referred to as wetted parts), which are widely accepted by life scientists, as they are known for optimum inertness to biological samples, and ensure best compatibility to common samples and solvents over a wide pH range. Explicitly, the complete flow path is free from stainless steel and free from other alloys containing metals such as iron, nickel, cobalt, chromium, molybdenum or copper, which can interfere with biological samples. The flow downstream of the sample introduction contains no metals whatsoever.

However, there are no materials that combine suitability for versatile HPLC instrumentation (valves, capillaries, springs, pump heads, flow cells etc.) with complete compatibility with all possible chemicals and application conditions. This section recommends the preferred solvents. Chemicals that are known to cause issues should be avoided, or exposure should be minimized, for example, for short-term cleaning procedures. After potentially aggressive chemicals have been used, the system should be flushed with compatible standard HPLC solvents.

PEEK

PEEK (Polyether-Ether Ketones) combines excellent properties with regard to biocompatibility, chemical resistance, mechanical and thermal stability and is therefore the material of choice for biochemical instrumentation. It is stable in the specified pH range, and inert to many common solvents. There is still a number of known incompatibilities with chemicals such as chloroform, methylene chloride, THF, DMSO, strong acids (nitric acid > 10 %, sulphuric acid > 10 %, sulfonic acids, trichloroacetic acid), halogenes or aqueous halogene solutions, phenol and derivatives (cresols, salicylic acid etc.).

When used above room temperature, PEEK is sensitive to bases and various organic solvents, which can cause it to swell. As normal PEEK capillaries are very sensitive to high pressure, especially under such conditions, Agilent uses stainless-steel clad PEEK capillaries to keep the flow path free of steel and to ensure pressure stability to at least 600 bar. If in doubt, consult the available literature about the chemical compatibility of PEEK.

Titanium

Titanium is highly resistant to oxidizing acids (for example, nitric, perchloric and hypochlorous acid) over a wide range of concentrations and temperatures. This is due to a thin oxide layer on the surface, which is stabilized by oxidizing compounds. Reducing acids (for example, hydrochloric, sulfuric and phosphoric acid) can cause slight corrosion, which increases with acid concentration and temperature. For example, the corrosion rate with 3 % HCl (about pH 0.1) at room temperature is about 13 µm/year. At room temperature, titanium is resistant to concentrations of about 5 % sulfuric acid (about pH 0.3). The addition of nitric acid to hydrochloric or sulfuric acids significantly reduces corrosion rates. Titanium is subject to corrosion in anhydrous methanol, which can be avoided by adding a small amount of water (about 3 %). Slight corrosion is possible with ammonia > 10 %.

Fused silica

Fused silica is inert against all common solvents and acids except hydrofluoric acid. It is corroded by strong bases and should not be used above pH 12 at room temperature. The corrosion of flow cell windows can negatively affect measurement results. For a pH greater than 12, the use of flow cells with sapphire windows is recommended.

Gold

Gold is inert to all common HPLC solvents, acids and bases within the specified pH range. It can be corroded by complexing cyanides and concentrated acids like aqua regia (a mixture of concentrated hydrochloric and nitric acid).

Zirconium Oxide

Zirconium Oxide (ZrO_2) is inert to almost all common acids, bases and solvents. There are no documented incompatibilities for HPLC applications.

Platinum/Iridium

Platinum/Iridium is inert to almost all common acids, bases and solvents. There are no documented incompatibilities for HPLC applications.

PTFE

PTFE (polytetrafluorethen) is inert to almost all common acids, bases and solvents. There are no documented incompatibilities for HPLC applications.

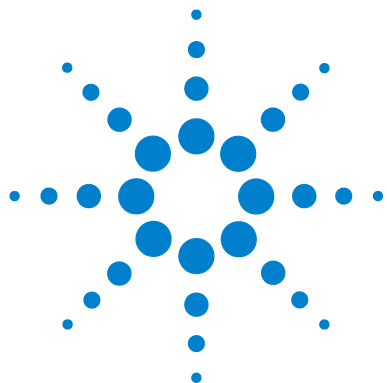
Sapphire, Ruby and Al_2O_3 -based ceramics

Sapphire, ruby and ceramics based on Al_2O_3 are inert to almost all common acids, bases and solvents. There are no documented incompatibilities for HPLC applications.

Data above were collected from external resources and are meant as a reference. Agilent cannot guarantee the completeness and correctness of such information. Information can also not be generalized due to catalytic effects of impurities like metal ions, complexing agents, oxygen etc. Most data available refers to room temperature (typically 20 – 25 °C, 68 – 77 °F). If corrosion is possible, it usually increases at higher temperatures. If in doubt, consult additional resources.

4 Using the Fraction Collector

Solvent Information



5

How to Optimize the Fraction Collector

Optimizing Fraction Collection [90](#)

This chapter provides information on how to optimize the module.



Optimizing Fraction Collection

Time-based fraction collection

- *Time slices* must have a length of at least 0.05 min.
- Set *# of Fractions* such that length of resulting fractions is at least 0.05 min.

Peak-based fraction collection

- Set threshold and slope values such that length of fractions is at least 0.05 min.
- Unresolved peaks can be separated using appropriate threshold and slope values. If two unresolved peaks are to be collected as one fraction, collect based on threshold only.
- If the baseline of the chromatogram is below or above 0 mAU, this offset is not accounted for when triggering peaks using a threshold value. The threshold value is always added to 0 mAU.



6 Troubleshooting and Diagnostics

Overview of the Module's Indicators and Test Functions [92](#)

Status Indicators [93](#)

Power Supply Indicator [93](#)

Module Status Indicator [94](#)

Agilent Lab Advisor Software [95](#)

Overview about the troubleshooting and diagnostic features.



Overview of the Module's Indicators and Test Functions

Status Indicators

The module is provided with two status indicators which indicate the operational state (prerun, run, and error states) of the module. The status indicators provide a quick visual check of the operation of the module.

Error Messages

In the event of an electronic, mechanical or hydraulic failure, the module generates an error message in the user interface. For each message, a short description of the failure, a list of probable causes of the problem, and a list of suggested actions to fix the problem are provided (see chapter Error Information).

Maintenance Functions

The maintenance functions position the transport unit and needle carrier assembly on certain positions for maintenance, homing, parking or delay calibration (see [“Maintenance Functions”](#) on page 118).

Transport Unit Self Alignment

The transport unit alignment with the transport unit and the well-plate tray is required to compensate for larger deviations in positioning the needle carrier assembly.

The transport unit self alignment is required after disassembling the system or when you exchange the transport unit, the needle carrier assembly or the MTP main board. This function is in the Instant Pilot and Lab Advisor.

Status Indicators

Two status indicators are located on the front of the module. The lower left indicates the power supply status, the upper right indicates the module status.

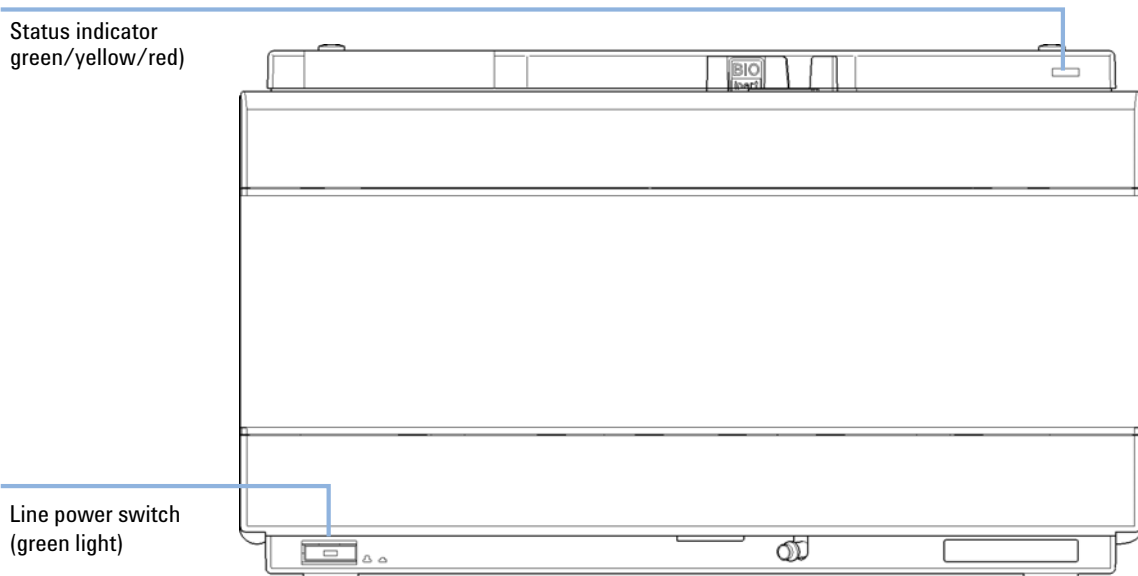


Figure 26 Location of status indicators

Power Supply Indicator

The power supply indicator is integrated into the main power switch. When the indicator is illuminated (*green*) the power is *ON*.

Module Status Indicator

The module status indicator indicates one of six possible module conditions:

- When the status indicator is *OFF* (and power switch light is on), the module is in a *prerun* condition, and is ready to begin an analysis.
- A *green* status indicator, indicates the module is performing an analysis (*run mode*).
- A *yellow* indicator indicates a *not-ready* condition. The module is in a not-ready state when it is waiting for a specific condition to be reached or completed (for example, immediately after changing a set point), or while a self-test procedure is running.
- An *error* condition is indicated when the status indicator is *red*. An error condition indicates the module has detected an internal problem which affects correct operation of the module. Usually, an error condition requires attention (e.g. leak, defective internal components). An error condition always interrupts the analysis.

If the error occurs during analysis, it is propagated within the LC system, i.e. a red LED may indicate a problem of a different module. Use the status display of your user interface for finding the root cause/module of the error.

- A *blinking* indicator indicates that the module is in resident mode (e.g. during update of main firmware).
- A *fast blinking* indicator indicates that the module is in a low-level error mode. In such a case try to re-boot the module or try a cold-start (see [“Special Settings”](#) on page 205). Then try a firmware update (see [“Replacing Module Firmware”](#) on page 162). If this does not help, a main board replacement is required.

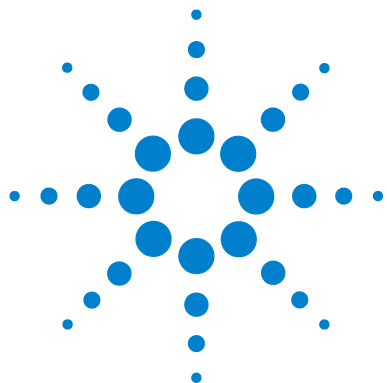
Agilent Lab Advisor Software

The Agilent Lab Advisor software is a standalone product that can be used with or without data system. Agilent Lab Advisor software helps to manage the lab for high quality chromatographic results and can monitor in real time a single Agilent LC or all the Agilent GCs and LCs configured on the lab intranet.

Agilent Lab Advisor software provides diagnostic capabilities for all Agilent 1200 Infinity Series modules. This includes diagnostic capabilities, calibration procedures and maintenance routines for all the maintenance routines.

The Agilent Lab Advisor software also allows users to monitor the status of their LC instruments. The Early Maintenance Feedback (EMF) feature helps to carry out preventive maintenance. In addition, users can generate a status report for each individual LC instrument. The tests and diagnostic features as provided by the Agilent Lab Advisor software may differ from the descriptions in this manual. For details refer to the Agilent Lab Advisor software help files.

The Instrument Utilities is a basic version of the Lab Advisor with limited functionality required for installation, use and maintenance. No advanced repair, troubleshooting and monitoring functionality is included.



7 Error Information

What Are Error Messages	99
General Error Messages	100
Timeout	100
Shutdown	101
Remote Timeout	101
Lost CAN Partner	102
Leak Sensor Short	102
Leak Sensor Open	103
Compensation Sensor Open	103
Compensation Sensor Short	104
Fan Failed	104
Leak	105
Open Cover	105
Module Error Messages	106
Exhaust Fan Failed	106
Front Door Error	106
Side Door Error	107
Arm Movement Failed or Arm Movement Timeout	107
Needle to Needle Rinse / Funnel Position Failed	108
Needle Carrier Failed	109
Missing Vial or Missing Well-plate	109
Calib delay vol two peaks	110
Valve Switch Failed	110
Adapter Required	111
Funnel Not Supported	111
Pusher Wrong or Defect	112
Wrong or Missing Needle (Analytical Scale)	112
Initialization Failed	112



Motor Temperature	113
Vessel Stuck to Needle	114
Cluster Partner Lost During Analysis	114
Movement to Next Position Failed	115
Could Not Find a Valid Next Position	115

This chapter describes the meaning of error messages, and provides information on probable causes and suggested actions how to recover from error conditions.

What Are Error Messages

Error messages are displayed in the user interface when an electronic, mechanical, or hydraulic (flow path) failure occurs which requires attention before the analysis can be continued (for example, repair, or exchange of consumables is necessary). In the event of such a failure, the red status indicator at the front of the module is switched on, and an entry is written into the module logbook.

General Error Messages

General error messages are generic to all Agilent series HPLC modules and may show up on other modules as well.

Timeout

Error ID: 0062

The timeout threshold was exceeded.

Probable cause

- 1** The analysis was completed successfully, and the timeout function switched off the module as requested.
- 2** A not-ready condition was present during a sequence or multiple-injection run for a period longer than the timeout threshold.

Suggested actions

Check the logbook for the occurrence and source of a not-ready condition. Restart the analysis where required.

Check the logbook for the occurrence and source of a not-ready condition. Restart the analysis where required.

Shutdown

Error ID: 0063

An external instrument has generated a shutdown signal on the remote line.

The module continually monitors the remote input connectors for status signals. A LOW signal input on pin 4 of the remote connector generates the error message.

Probable cause	Suggested actions
1 Leak detected in another module with a CAN connection to the system.	Fix the leak in the external instrument before restarting the module.
2 Leak detected in an external instrument with a remote connection to the system.	Fix the leak in the external instrument before restarting the module.
3 Shut-down in an external instrument with a remote connection to the system.	Check external instruments for a shut-down condition.

Remote Timeout

Error ID: 0070

A not-ready condition is still present on the remote input. When an analysis is started, the system expects all not-ready conditions (for example, a not-ready condition during detector balance) to switch to run conditions within one minute of starting the analysis. If a not-ready condition is still present on the remote line after one minute the error message is generated.

Probable cause	Suggested actions
1 Not-ready condition in one of the instruments connected to the remote line.	Ensure the instrument showing the not-ready condition is installed correctly, and is set up correctly for analysis.
2 Defective remote cable.	Exchange the remote cable.
3 Defective components in the instrument showing the not-ready condition.	Check the instrument for defects (refer to the instrument's documentation).

Lost CAN Partner

Error ID: 0071

During an analysis, the internal synchronization or communication between one or more of the modules in the system has failed.

The system processors continually monitor the system configuration. If one or more of the modules is no longer recognized as being connected to the system, the error message is generated.

Probable cause

- 1 CAN cable disconnected.
- 2 Defective CAN cable.
- 3 Defective main board in another module.

Suggested actions

- Ensure all the CAN cables are connected correctly.
 - Ensure all CAN cables are installed correctly.
- Exchange the CAN cable.
- Switch off the system. Restart the system, and determine which module or modules are not recognized by the system.

Leak Sensor Short

Error ID: 0082

The leak sensor in the module has failed (short circuit).

The current through the leak sensor is dependent on temperature. A leak is detected when solvent cools the leak sensor, causing the leak sensor current to change within defined limits. If the current increases above the upper limit, the error message is generated.

Probable cause

- 1 Defective leak sensor.
- 2 Leak sensor incorrectly routed, being pinched by a metal component.

Suggested actions

- Please contact your Agilent service representative.
- Please contact your Agilent service representative.

Leak Sensor Open

Error ID: 0083

The leak sensor in the module has failed (open circuit).

The current through the leak sensor is dependent on temperature. A leak is detected when solvent cools the leak sensor, causing the leak-sensor current to change within defined limits. If the current falls outside the lower limit, the error message is generated.

Probable cause	Suggested actions
1 Leak sensor not connected to the main board.	Please contact your Agilent service representative.
2 Defective leak sensor.	Please contact your Agilent service representative.
3 Leak sensor incorrectly routed, being pinched by a metal component.	Please contact your Agilent service representative.

Compensation Sensor Open

Error ID: 0081

The ambient-compensation sensor (NTC) on the main board in the module has failed (open circuit).

The resistance across the temperature compensation sensor (NTC) on the main board is dependent on ambient temperature. The change in resistance is used by the leak circuit to compensate for ambient temperature changes. If the resistance across the sensor increases above the upper limit, the error message is generated.

Probable cause	Suggested actions
1 Defective main board.	Please contact your Agilent service representative.

Compensation Sensor Short

Error ID: 0080

The ambient-compensation sensor (NTC) on the main board in the module has failed (short circuit).

The resistance across the temperature compensation sensor (NTC) on the main board is dependent on ambient temperature. The change in resistance is used by the leak circuit to compensate for ambient temperature changes. If the resistance across the sensor falls below the lower limit, the error message is generated.

Probable cause

- 1 Defective main board.

Suggested actions

Please contact your Agilent service representative.

Fan Failed

Error ID: 0068

The cooling fan in the module has failed.

The hall sensor on the fan shaft is used by the main board to monitor the fan speed. If the fan speed falls below a certain limit for a certain length of time, the error message is generated.

Depending on the module, assemblies (e.g. the lamp in the detector) are turned off to assure that the module does not overheat inside.

Probable cause

- 1 Fan cable disconnected.
- 2 Defective fan.
- 3 Defective main board.

Suggested actions

Please contact your Agilent service representative.
Please contact your Agilent service representative.
Please contact your Agilent service representative.

Leak

Error ID: 0064

A leak was detected in the module.

The signals from the two temperature sensors (leak sensor and board-mounted temperature-compensation sensor) are used by the leak algorithm to determine whether a leak is present. When a leak occurs, the leak sensor is cooled by the solvent. This changes the resistance of the leak sensor which is sensed by the leak-sensor circuit on the main board.

Probable cause

- 1 Loose fittings.
- 2 Broken capillary.

Suggested actions

- Ensure all fittings are tight.
- Exchange defective capillaries.

Open Cover

Error ID: 0205

The top foam has been removed.

Probable cause

- 1 Foam not activating the sensor.
- 2 Dirty or defective sensor.

Suggested actions

- Please contact your Agilent service representative.
- Please contact your Agilent service representative.

Module Error Messages

Exhaust Fan Failed

Error ID: 4456, 4457

The exhaust fan in the module has failed.

The hall sensor on the fan shaft is used by the main board to monitor the fan speed. If the fan speed falls below a certain value the error message is generated and the module shuts down.

Probable cause	Suggested actions
1 Fan cable disconnected.	Please contact your Agilent service representative.
2 Defective fan.	Please contact your Agilent service representative.
3 Defective main board.	Please contact your Agilent service representative.

Front Door Error

Error ID: 4750

The front door and/or the SLF board are damaged.

Probable cause	Suggested actions
1 The flat ribbon cable from MTP main board to the SLF board is not connected correctly.	Please contact your Agilent service representative.
2 The sensor on the SLF board is defective.	Please contact your Agilent service representative.
3 The door is bent or the magnet is misplaced/broken.	Change the side door.
4 Defective MTP main board.	Please contact your Agilent service representative.

Side Door Error

Error ID: 4750

The side door and/or the MTP board are damaged. This error message is not displayed before the initialization is finished.

Probable cause	Suggested actions
1 The side door is not installed.	Install the side door.
2 The door is bent or the magnet is misplaced/broken.	Change the side door.
3 The sensor on the MTP board is defective.	Please contact your Agilent service representative.
4 Excessive weight on top of the fraction collector (see also " Initialization Failed " on page 112)	Check stack configuration and reduce weight on top of the fraction collector.

Arm Movement Failed or Arm Movement Timeout

Error ID: 4002

The transport assembly was unable to complete a movement in one of the axes.

The processor defines a certain time window for the successful completion of a movement in any particular axis. The movement and position of the transport assembly is monitored by the encoders on the stepper motors. If the processor does not receive the correct position information from the encoders within the time window, the error message is generated.

Axes identification:

- Arm Movement 0 Failed: X-axis.
- Arm Movement 1 Failed: Z-axis.
- Arm Movement 2 Failed: Theta (needle carrier rotation).

Probable cause	Suggested actions
1 Mechanical obstruction.	Ensure unobstructed movement of the transport assembly.
2 High friction in the transport assembly.	Please contact your Agilent service representative.
3 Defective motor assembly.	Please contact your Agilent service representative.
4 Defective sample transport assembly flex board.	Please contact your Agilent service representative.
5 Defective main board.	Please contact your Agilent service representative.

Needle to Needle Rinse / Funnel Position Failed

Error ID: 4955, 4980, 4981-4990

The needle failed to reach the needle rinse / funnel position.

The position of the needle is monitored by a position encoder on the needle carrier. If the needle fails to reach the end point, or if the encoder fails to recognize the needle carrier movement, the error message is generated.

Probable cause	Suggested actions
1 Bad sample transport unit alignment	Do a self-alignment
2 Bent needle.	Check and exchange the needle assembly if necessary.
3 Missing needle.	Check and exchange the needle assembly if necessary.
4 Blocked rinse or funnel position.	Clean or change the funnel or rinse port assembly if necessary.
5 Defective needle carrier assembly.	Exchange the needle carrier assembly.
6 Disconnected needle carrier connector.	Connect needle carrier connector correctly.
7 Defective MTP main board.	Please contact your Agilent service representative.

Needle Carrier Failed

Error ID:

The needle carrier on the transport unit assembly failed to move correctly.

Probable cause	Suggested actions
1 Defective position sensor in the needle carrier assembly.	Exchange the needle carrier assembly.
2 Bad needle carrier positioning in X or Theta.	Perform a self-alignment.
3 Disconnected needle carrier connector.	Connect needle carrier connector correctly.
4 Defective MTP main board.	Please contact your Agilent service representative.
5 Defective Z-motor.	Please contact your Agilent service representative.

Missing Vial or Missing Well-plate

Error ID:

No vial or well-plate was found in the position defined in the method or sequence.

When the needle carrier moves to a vial or well-plate and the needle is lowered into the vial or well-plate, the position of the needle is monitored by an encoder behind the vial pusher. If no vial or well-plate is present, the encoder detects an error and the message “missing vial or well plate” is generated.

Probable cause	Suggested actions
1 No vial in the position defined in the method or sequence.	Install the sample vial in the correct position, or edit the method or sequence accordingly.
2 Defective needle carrier assembly.	Exchange the needle carrier assembly.
3 Defective transport unit assembly flex board.	Please contact your Agilent service representative.
4 Defective MTP main board.	Please contact your Agilent service representative.

Calib delay vol two peaks

Error ID: 4759

Two peaks have been detected during the delay calibration.

Probable cause	Suggested actions
1 Wrong sample has been used for the delay calibration.	Check method and delay calibration procedure.
2 Wrong method has been used for the delay calibration.	Check method and delay calibration procedure.
3 Air bubbles are in the flow path.	Check flow path for leaks and air bubbles.

Valve Switch Failed

Error ID: 4959

If multiple fraction collectors are configured, an external valve G1160A is used to switch between the fraction collectors.

The error message is displayed if the external valve failed to switch to next position.

Probable cause	Suggested actions
1 Valve is blocked. Possible if eluents with highly concentrated electrolytes are used.	Purge valve to dissolve crystals.
2 Power cord for the valve is not connected.	Check power cord connection.
3 Valve drive or valve electronics are defective.	• Synchronize the valve. • Exchange the valve.

Adapter Required

Error ID: 4961

The wellplate adapter is required for the following operation.

Probable cause	Suggested actions
1 Wellplate adapter not attached.	Attach wellplate adapter.
2 Cable between needle carrier assembly to transport unit assembly disconnected or defective.	Check cable between needle carrier assembly to transport unit assembly.
3 Needle carrier assembly defective.	Exchange needle carrier assembly.

Funnel Not Supported

Error ID: 4962

Funnels are only supported for the analytical scale fraction collector. Consequently the error message is displayed, if a tray with funnels configured for the preparative or the micro collector/spotter.

Probable cause	Suggested actions
1 Fraction collector type and tray not compatible.	- Check your tray configuration. Select Instrument (Agilent 1100/1200 System) > More Agilent 1100/1200 Fraction Collector > Configuration. Funnel trays are only supported for the analytical scale fraction collector. - Check your 1260 Infinity System Configuration. Select Instrument (Agilent 1100/1200 System) > More Agilent 1100/1200 Fraction Collector > Configure 1100/1200 Access... and verify, that an analytical scale fraction collector is configured.

Pusher Wrong or Defect

Error ID: 4965

Probable cause

- 1 Wellplate adapter not attached.
- 2 Needle carrier assembly defective.

Suggested actions

- Attach wellplate adapter.
- Exchange needle carrier assembly.

Wrong or Missing Needle (Analytical Scale)

Error ID: 4966

Probable cause

- 1 No needle installed.
- 2 Short needle for high flow rates installed, but the wellplate adapter hasn't been removed.

Suggested actions

- Check which needle has been installed.
- Remove wellplate adapter if short needle for semi-preparative operation is installed.

Initialization Failed

Error ID: 4950

The fraction collector failed to complete initialization correctly. The fraction collector initialization procedure moves the needle arm and transport assembly to their home positions in a predefined routine. During initialization, the processor monitors the position sensors and motor encoders to check for correct movement. If one or more of the movements is not successful, or is not detected, the error message is generated.

Probable cause

- 1 Transport unit not aligned correctly
- 2 Mechanical obstruction.
- 3 Defective transport assembly flex board.

Suggested actions

- Perform an auto-alignment.
- Ensure unobstructed movement of the transport assembly.
- Please contact your Agilent service representative.

Probable cause	Suggested actions
4 Defective MTP main board.	Please contact your Agilent service representative.
5 Excessive weight on top of the fraction collector (see also “ Side Door Error ” on page 107)	Check stack configuration and reduce weight on top of the fraction collector.

Motor Temperature

Error ID: 4027, 4040, 4261, 4451

One of the motors of the transport assembly has drawn excessive current, causing the motor to become too hot. The processor has switched off the motor to prevent damage to the motor.

Motor identification:

- Motor 0 temperature: X-axis motor.
- Motor 1 temperature: Z-axis motor.
- Motor 2 temperature: Theta motor.

The processor monitors the current drawn by each motor and the time the motor is drawing current. The current drawn by the motors is dependent on the load on each motor (friction, mass of components etc.). If the current drawn is too high, or the time the motor draws current is too long, the error message is generated.

Probable cause	Suggested actions
1 Mechanical obstruction.	Ensure unobstructed movement of the transport assembly.
2 High friction in the transport assembly.	Please contact your Agilent service representative.
3 Motor belt tension too high.	Switch off the module at the power switch. Wait at least 10 minutes before switching on again.
4 Defective motor.	Please contact your Agilent service representative.
5 Defective transport assembly flex board.	Please contact your Agilent service representative.

Vessel Stuck to Needle

Error ID: 4453

The vessel sticks to the needle when the needle moves up.

Probable cause

- 1** Closing mat to rigid/thick.
- 2** Bad X or Theta positioning and the needle sticks into the wall between two holes.
- 3** Defective encoder on the needle carrier assembly.

Suggested actions

- Check that the closing mat is not too thick.
- Please contact your Agilent service representative.
- Please contact your Agilent service representative.

Cluster Partner Lost During Analysis

Error ID:

There was a problem with the inter module communication.

Probable cause

- 1** Disconnected or defective CAN cable.
- 2** Disconnected or defective 24V-CAN-DC-OUT cable for an external valve.
- 3** Defective UIB, external Valve or MTP board.

Suggested actions

- Check the interconnection between the modules.
- Reconnect the UIB / Valve. Start a test analysis/run.
- Switch power off / on (complete system off, then on). Start a test analysis / run.
 - Please contact your Agilent service representative.

Movement to Next Position Failed

Error ID: 4957

The transport mechanism detected an unexpected situation during the movement to the next fraction position.

Probable cause

- 1 Mismatch between tray configuration and the loading of the trays, e.g. 4 well plates are configured in the UI, but only three are loaded, or shallow plates are configured, but deep well-plates are used, or tube height doesn't match the configuration.

Suggested actions

Check the configuration and the loading of the tray.

Could Not Find a Valid Next Position

Error ID: 4958

There has been more fractions than fraction positions.

Probable cause

- 1 Unexpected number of fractions.
- 2 Start of an analysis / sequence / run without changing of the tray of the previous run.

Suggested actions

- If possible use a tray with more positions.
 - Add an additional G5664 to the system.
- Ensure that there are enough fraction positions for the complete analysis / sequence / run.

7 Error Information

Module Error Messages



8 Test Functions

Maintenance Functions [118](#)

Transport Unit Self Alignment [120](#)

Step Commands [122](#)

Troubleshooting [123](#)

This chapter describes the module's built in test functions.



Maintenance Functions

Some maintenance procedures require the needle arm and needle carrier to be moved to specific positions to enable easy access to components. The maintenance functions move these assemblies into the appropriate maintenance position. In the ChemStation, the **Fraction Collector Maintenance Positions** can be selected from the **Maintenance** menu in the **Diagnosis** view (see [Figure 27](#) on page 118). In the Instant Pilot, the functions can be selected in the **Test** screens of the fraction collector.

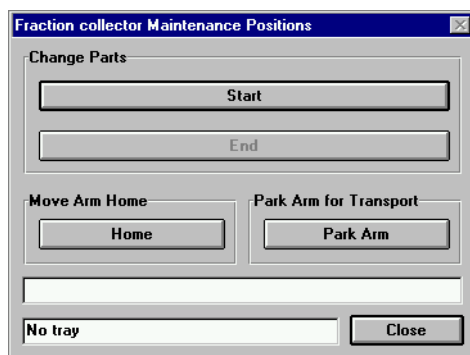


Figure 27 Fraction Collector Maintenance Positions dialog box

Change Parts

If you click **Start** the transport unit will move upwards, the needle carrier assembly will move to the front center and then turn off the theta motor to allow free rotation of the theta arm. This position enables easy access to the transport unit to:

- change the needle assembly
- change the needle/capillary carrier assembly,
- change the valve to needle tubing
- change the diverter valve

After the maintenance or repair task has been finished, click **End** to move the transport assembly to the **Home** position.

Home Position

This maintenance function moves the arm up and to the right rear for better access and exchange of the trays.

Park Arm

This maintenance position moves the arm to the park position at the upper rear left side of the tray for transporting or shipping the fraction collector.

Transport Unit Self Alignment

The transport unit alignment is required to compensate for larger deviations in positioning the needle carrier assembly. This might be necessary after disassembling the system or when you exchange the transport unit, the needle carrier assembly or the MTP main board. This function is in the **Diagnose** screen of the ChemStation, the Instant Pilot and Lab Advisor.

When

- After disassembling the module
- When exchanging:
 - the transport unit
 - the needle/capillary carrier assembly
 - the MTP main board

CAUTION

Damage to the module

If the Transport Unit Self Alignment is started with well plates on the tray, the alignment procedure is aborted WITHOUT error message.

- The sample transport self alignment requires the 4-well-plate tray (p/n G1364-84521).
- Remove all wellplates.

- 1 If the transport unit has been exchanged or if it is strongly misaligned, set the 8-bit configuration switch to the Forced Cold Start Configuration, see [Table 26](#) on page 205.
- 2 Install 4-well-plate tray (p/n G1364-84521).

NOTE

Remove all plates!

- 3 Ensure that the wellplate adapter is correctly assembled.
- 4 Select the **Maintenance** menu in the **Diagnosis** view of the Agilent ChemStation.

- 5 In the menu choose **Fraction Collector > Transport Alignment...** to start the automated procedure. The Transport Alignment Procedure takes approximately 10-15 minutes.
- 6 Set the 8-bit configuration switch to the default setting, see [Table 20](#) on page 203 .

NOTE

If the Transport Unit Self Alignment is started with well plates on the tray, the alignment procedure is aborted without any error message.

Step Commands

Some movements of the fraction collection sequence can be done under manual control. This is useful during troubleshooting where close observation of each of the fraction collection step is required to confirm a specific failure mode or verify successful completion of a repair.

Each step command actually consists of a series of individual commands which move the fraction collector components to predefined positions enabling the specific step to be done.

In Lab Advisor the step commands can be selected from the Tools screen.

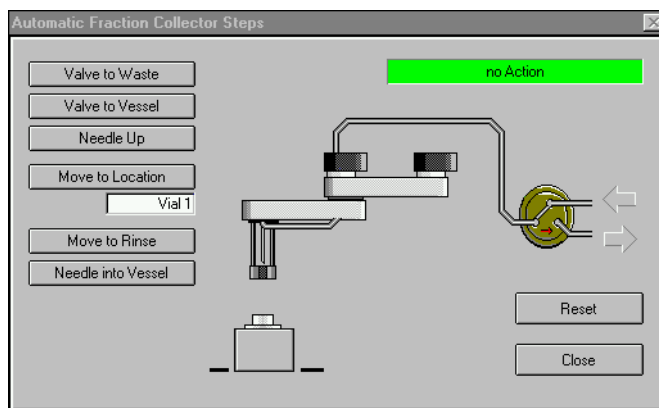


Figure 28 Fraction Collector Step Commands

Table 15 Step Commands

Step	Action	Comments
Needle Up	Lifts the needle arm to the upper position.	Command also switches the diverter valve to waste if it is not already in that position.
Needle into vessel	Lowest the needle into the specified vessel.	
Needle to rinse / flush port	Moves the needle to the rinse / flush port.	
Switch valve to needle	Switches the diverter valve to the needle.	
Switch valve to waste	Switches the diverter valve from needle to waste.	

Troubleshooting

If the fraction collector is unable to perform a specific step due to a hardware failure, an error message is generated. You can use the step commands to perform a fraction collection sequence, and observe how the fraction collector responds to each command.



9 Maintenance

Introduction to Maintenance	126
Cautions and Warnings	127
Cleaning the Module	129
Overview of Maintenance	130
Replacing the Inlet/Waste Tubings	132
Replacing the Valve to Needle Tubing	136
Exchanging the Needle Assembly	140
Exchanging the Needle/Capillary Carrier Assembly	143
Exchanging the Diverter Valve	148
Exchanging the Internal Tray	152
Repairing or Exchanging a Funnel of the Internal Tray	154
Exchanging the Leak Sensor	157
Exchanging or Installing the (Optional) Interface Board	160
Replacing Module Firmware	162

This chapter describes the maintenance of the module.



Introduction to Maintenance

The module is designed for easy maintenance. Maintenance can be done from the front with module in place in the system stack.

NOTE

There are no serviceable parts inside.

Do not open the module.

Cautions and Warnings

WARNING

Module is partially energized when switched off, as long as the power cord is plugged in.

Risk of stroke and other personal injury. Repair work at the module can lead to personal injuries, e. g. shock hazard, when the module cover is opened and the instrument is connected to power.

- Never perform any adjustment, maintenance or repair of the module with the top cover removed and with the power cord plugged in.
 - The security lever at the power input socket prevents that the module cover is taken off when line power is still connected. Never plug the power line back in when cover is removed.
-

WARNING

Sharp metal edges

Sharp-edged parts of the equipment may cause injuries.

- To prevent personal injury, be careful when getting in contact with sharp metal areas.
-

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.
 - The volume of substances should be reduced to the minimum required for the analysis.
 - Do not operate the instrument in an explosive atmosphere.
-

CAUTION

Electronic boards and components are sensitive to electrostatic discharge (ESD). ESD can damage electronic boards and components.

- Be sure to hold the board by the edges, and do not touch the electrical components. Always use ESD protection (for example, an ESD wrist strap) when handling electronic boards and components.
-

CAUTION

Safety standards for external equipment

- If you connect external equipment to the instrument, make sure that you only use accessory units tested and approved according to the safety standards appropriate for the type of external equipment.
-

CAUTION

Sample degradation and contamination of the instrument

Metal parts in the flow path can interact with the bio-molecules in the sample leading to sample degradation and contamination.

- For bio-inert applications, always use dedicated bio-inert parts, which can be identified by the bio-inert symbol or other markers described in this manual.
 - Do not mix bio-inert and non-inert modules or parts in a bio-inert system.
-

NOTE

The electronics of the module will not allow operation of the module when the top cover and the top foam are removed. A safety light switch on the main board will inhibit the operation of the fan immediately. Voltages for the other electronic components will be turned off after 30 seconds. The status lamp will light up red and an error will be logged into the logbook of the user interface. Always operate the module with the top covers in place.

Cleaning the Module

To keep the module case clean, use a soft cloth slightly dampened with water, or a solution of water and mild detergent.

WARNING

Liquid dripping into the electronic compartment of your module can cause shock hazard and damage the module

- Do not use an excessively damp cloth during cleaning.
 - Drain all solvent lines before opening any connections in the flow path.
-

Overview of Maintenance

The procedures described in this section can be done with the fraction collector in place in the stack. These procedures can be done on a more frequent basis.

Table 16 Simple Repair Procedures

Procedure	Typical Frequency	Notes
Replacing the inlet / waste tubings	When worn out, when showing visual signs of damage, typically once per year.	See “Replacing the Inlet/Waste Tubings” on page 132
Replacing the valve to needle tubings	When worn out, when showing visual signs of damage, typically once per year	See “Replacing the Valve to Needle Tubing” on page 136
Exchanging the needle assembly	When needle shows indication of damage or blockage Or when using the short needle assembly for operation with high test tubes (45 mm>)	See “Exchanging the Needle Assembly” on page 140
Exchanging the needle/capillary carrier assembly	When the needle carrier is defective	See “Exchanging the Needle/Capillary Carrier Assembly” on page 143
Exchanging the diverter valve	When defective (internal / external leak, valve not switching any more)	See “Exchanging the Diverter Valve” on page 148
Exchanging the internal tray	When flow delay sensor defective	See “Exchanging the Internal Tray” on page 152
Repairing or exchanging a funnel of the internal tray or funnel tray	When defective (leaky, blocked or contaminated)	See “Repairing or Exchanging a Funnel of the Internal Tray” on page 154

Table 16 Simple Repair Procedures

Procedure	Typical Frequency	Notes
Exchanging the leak sensor	When defective	See “ Exchanging the Leak Sensor ” on page 157
Exchanging or installing the (optional) BCD board	When defective or new board installed	See “ Exchanging or Installing the (Optional) Interface Board ” on page 160

Replacing the Inlet/Waste Tubings

When	• When contaminated, worn out or visibly damaged • Typically once every year					
Parts required	<table><tr><th>p/n</th><th>Description</th></tr><tr><td>G5664-68712</td><td>Analytical tubing kit 0.25 mm i.d. PTFE-ESD</td></tr></table>	p/n	Description	G5664-68712	Analytical tubing kit 0.25 mm i.d. PTFE-ESD	
p/n	Description					
G5664-68712	Analytical tubing kit 0.25 mm i.d. PTFE-ESD					
Preparations	• Position the transport unit of the fraction collector in the Home position (see “Maintenance Functions” on page 118). • Remove all installed trays from the tray base. • Position the transport unit of the fraction collector in the Change Parts position (see “Maintenance Functions” on page 118). • Turn OFF the instrument. • Remove the rear end of the fraction collector’s waste tubing from the waste container, unscrew the front end of the fraction collector’s inlet tubing from the flow cell of the detector.					

WARNING

Personal injury

There is a risk of personal injury because the needle arm can move during fraction collector operation.

→ To avoid personal injury, keep fingers away from the needle area.

CAUTION

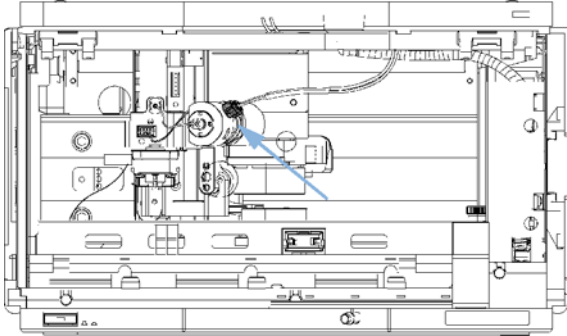
Liquid spills or fraction losses

Worn or damaged tubings can cause potential spills or lead to fraction losses.

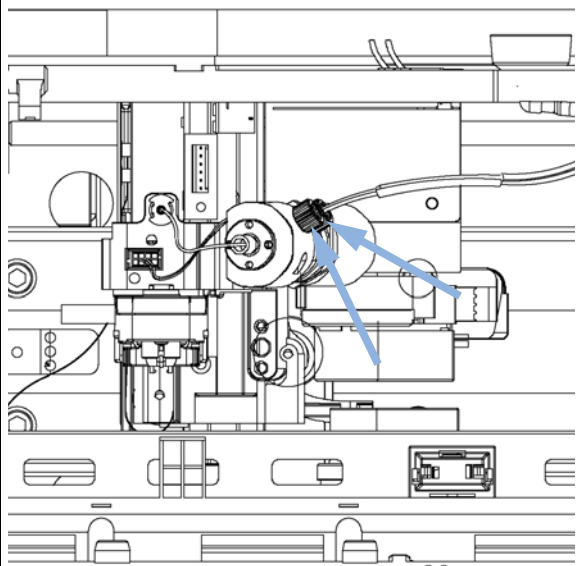
→ Explicitly follow the described installation procedures to maximize the lifetime of inlet / waste tubing assembly and the valve to needle tubing, and to avoid potential spills or fraction losses.

→ Regularly inspect the tubings and exchange them if they are worn out or show visible signs of damage.

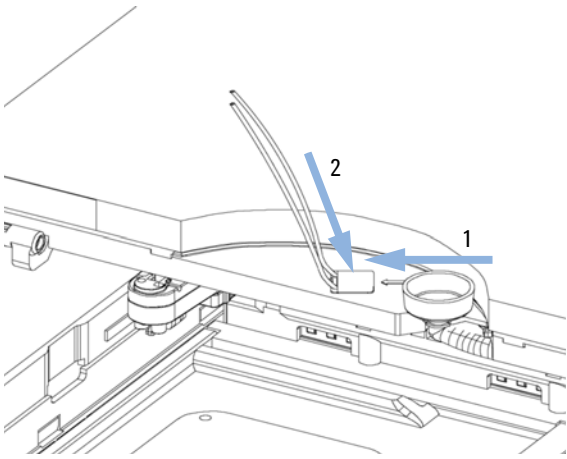
- 1** Locate the diverter valve with the finger tight fittings of the inlet / waste tubing assembly (the figure shows the open fraction collector seen from the front).



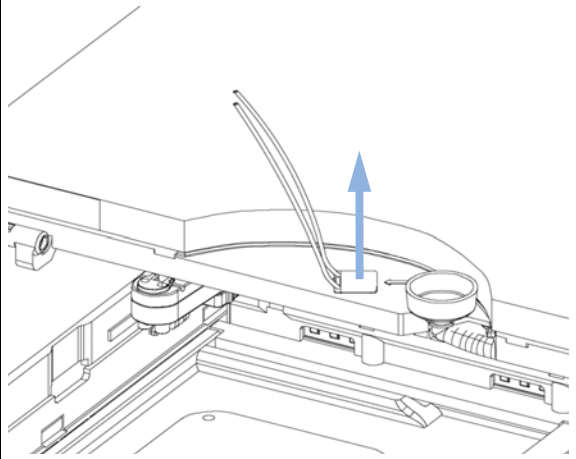
- 2** Unscrew the two finger-tight fittings of the inlet / waste tubing assembly at the diverter valve.



- 3** Unplug the inlet / waste tubing assembly from the bolt carrier (located on the top center of the front cover) by moving the snapper as indicated by the arrows. Slide the tubings out to the bottom.



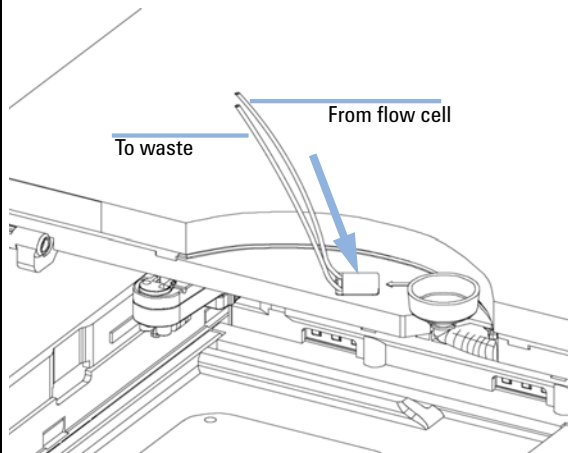
- 4** Install the new inlet / waste tubing assembly into the bolt carrier as shown below. Slide in the long ends of the tubings from bottom to top and let the snapper click into position.



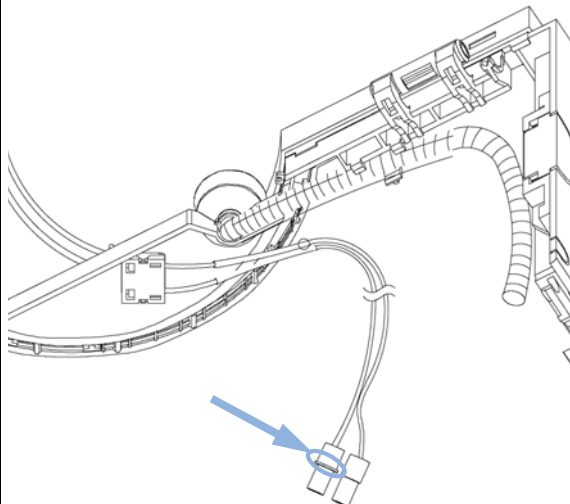
9 Maintenance

Replacing the Inlet/Waste Tubings

- 5** The rear tubing (with the label with the arrow) must be connected to the flow cell of the detector. The front tubing must be inserted into the waste container.



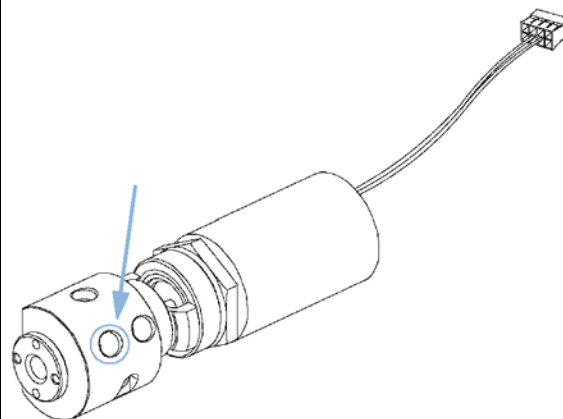
- 6** Connect the finger-tight fittings of the inlet / waste tubing assembly to the ports of the diverter valve. (View from the bottom)



NOTE

The tubings must not be bent up- or downwards. The cables must not be twisted.

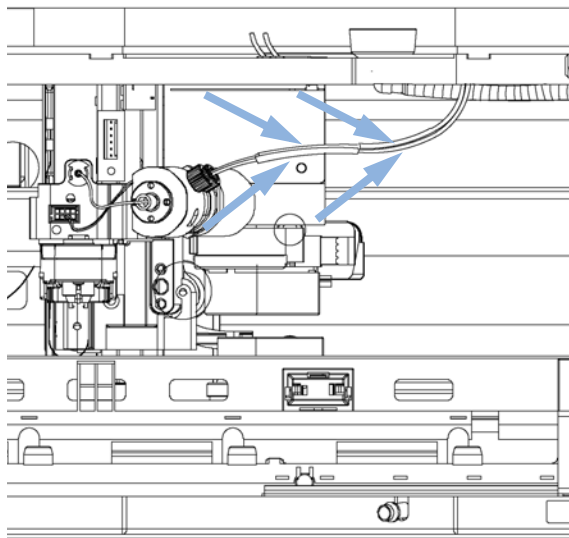
- 7** A color coded ring on one of the tubings and the valve body indicates, which cable belongs to which port.



NOTE

It is absolutely vital to connect these tubings as described, in order to maximize their lifetime and operating security.

8 IMPORTANT: The cables must run into the ports of the diverter valve in lines parallel to the horizon.



Next Steps:

- 9** Re-install the tray(s) in the tray base.
- 10** Start the instrument.
- 11** Close the front cover.

Replacing the Valve to Needle Tubing

When	• When contaminated, worn out or visibly damaged • Typically once every year	
Tools required	p/n	Description
	8710-1534	Wrench, open end, 4 mm
	8710-0510	Wrench open 1/4 — 5/16 inch
Parts required	p/n	Description
	G5664-68712	Analytical tubing kit 0.25 mm i.d. PTFE-ESD
Preparations	• Position the transport unit of the fraction collector in the Home position (see “Maintenance Functions” on page 118). • Remove all installed trays from the tray base. • Position the transport unit of the fraction collector in the Change Parts position (see “Maintenance Functions” on page 118) and turn off the instrument. • It might be more convenient to remove the needle from its carrier before unscrewing the needle tubing.	

WARNING

Personal injury

There is a risk of personal injury because the needle arm can move during fraction collector operation.

→ To avoid personal injury, keep fingers away from the needle area.

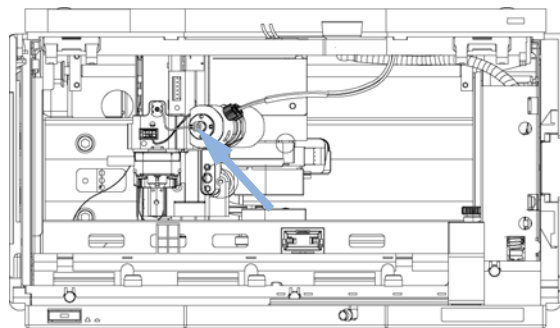
CAUTION

Liquid spills or fraction losses

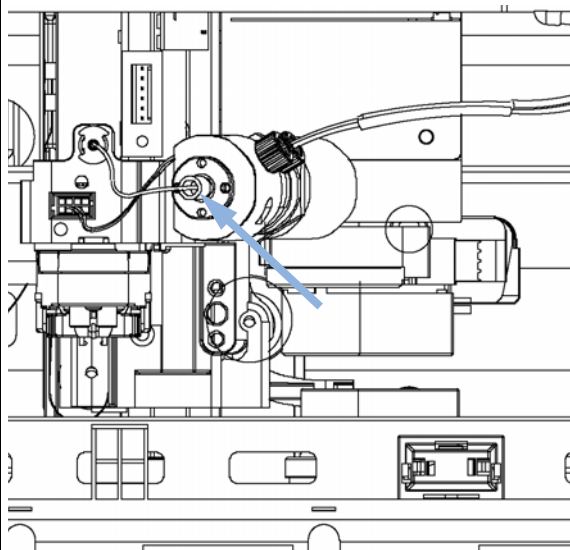
Worn or damaged tubings can cause potential spills or lead to fraction losses.

- Explicitly follow the described installation procedures to maximize the lifetime of inlet / waste tubing assembly and the valve to needle tubing, and to avoid potential spills or fraction losses.
- Regularly inspect the tubings and exchange them if they are worn out or show visible signs of damage.

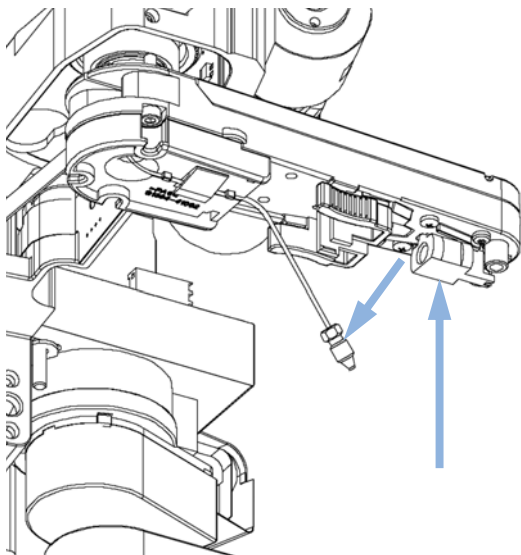
- 1** Locate the diverter valve with the finger-tight fittings of the valve to needle tubing assembly (the figure shows the open fraction collector seen from the front).



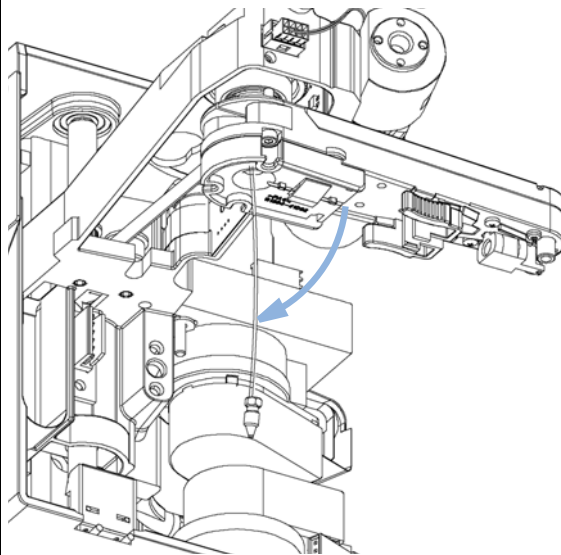
- 2** Unscrew the finger-tight fitting of the valve to needle tubing assembly at the diverter valve.



- 3** Using the 4 mm wrench and the 5/16" wrench for counter-holding unscrew the valve to needle tubing from the needle. (Viewed from the bottom)



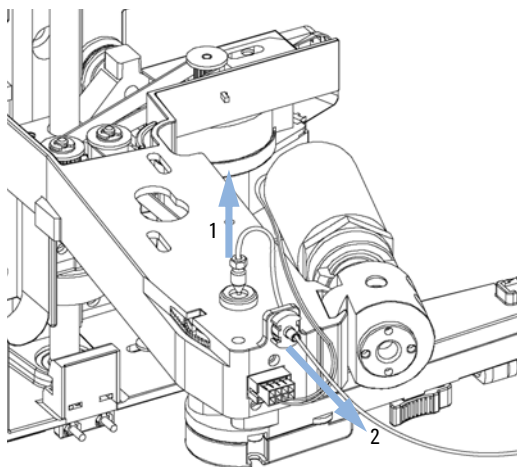
- 4** Un-clip the tubing from the bottom of the needle carrier assembly.



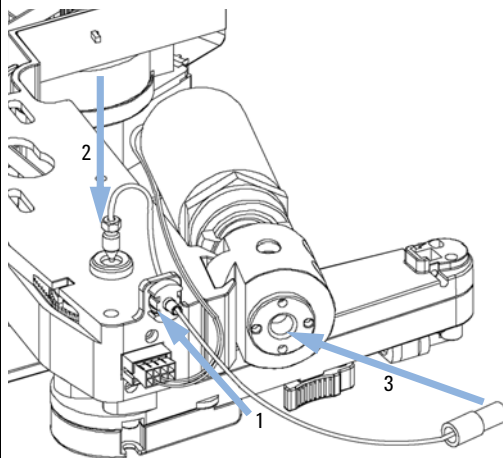
9 Maintenance

Replacing the Valve to Needle Tubing

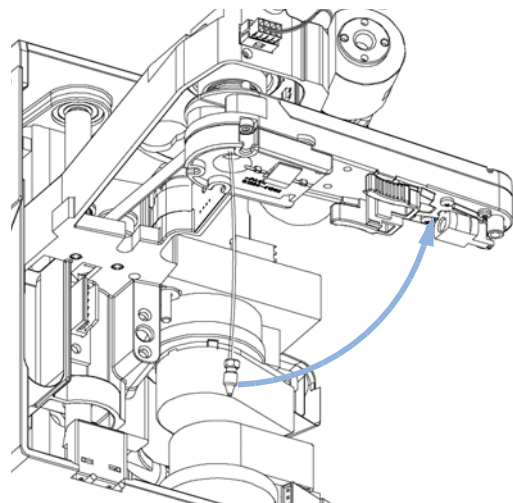
- 5** Slide the tubing through the hole in the needle carrier assembly (from bottom to top) and out of the holder in the z-arm assembly.



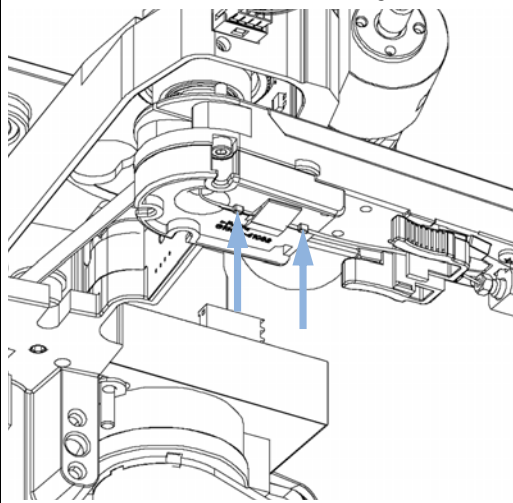
- 6** Install the new valve to needle tubing assembly by clipping it in to the holder in the z-arm assembly (1), Important!!! and slide it through the hole in the z-arm (2) and out on the bottom of the needle carrier assembly (top to bottom). Screw the finger-tight fitting into the port of the diverter valve (3).



- 7** Using the 4 mm wrench and the 5/16" wrench for counter-holding connect the valve to needle tubing to the needle. (Viewed from the bottom)



- 8** **IMPORTANT:** After fixing the screw clip the tubing into the guide on the bottom of the needle carrier assembly. It is absolutely vital that the tubing is installed as described, to maximize the lifetime of the tubing.



Next Steps:

- 9** Re-install the needle to the needle carrier assembly, if you previously removed it. Make sure to slide the needle all the way to the front of the needle carrier assembly (clicks into position).
- 10** Re-install the tray(s) in the tray base.
- 11** Start the instrument.
- 12** Close the front cover.

Exchanging the Needle Assembly

When	• When the needle is visibly damaged • When the needle is blocked or contaminated	
Tools required	p/n	Description
	8710-1534	Wrench, open end, 4 mm
	8710-0510	Wrench open 1/4 — 5/16 inch
Parts required	p/n	Description
	G5667-87200	Needle assembly (bio-inert)
Preparations	• Position the transport unit of the fraction collector in the Home position (see “Maintenance Functions” on page 118). • Remove all installed trays from the tray base. • Position the transport unit of the fraction collector in the Change Parts position (see “Maintenance Functions” on page 118) and turn off the instrument. • It might be more convenient to remove the needle from its carrier before unscrewing the needle tubing.	

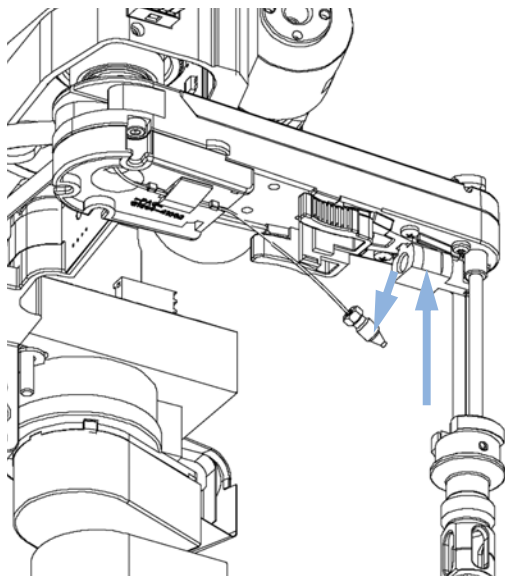
WARNING

Personal injury

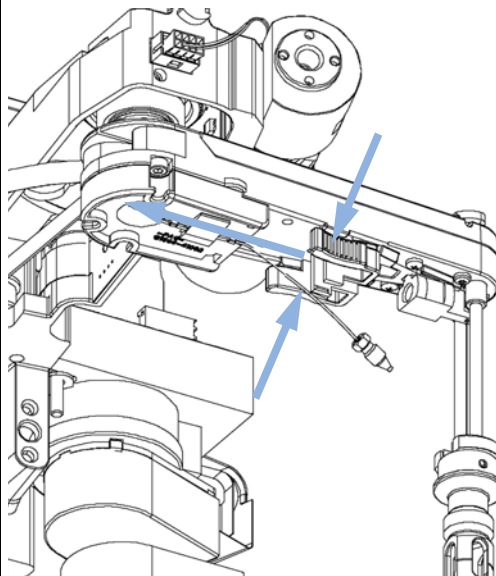
There is a risk of personal injury because the needle arm can move during fraction collector operation.

→ To avoid personal injury, keep fingers away from the needle area.

- 1** Use the 4 mm wrench and the 5/16" wrench for counter-holding to unscrew the valve-to-needle tubing from the needle assembly. (Viewed from the bottom).



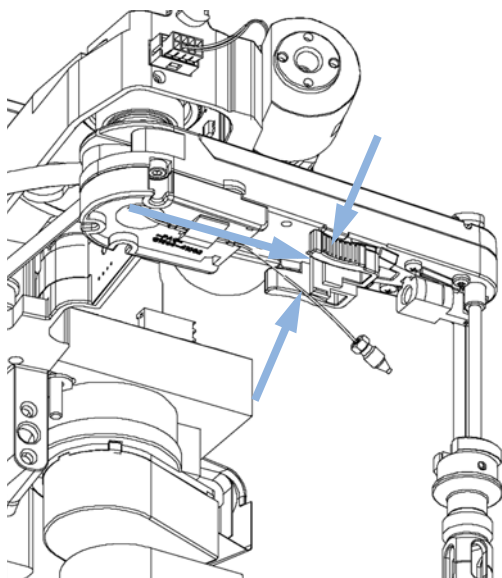
- 2** Remove the wellplate adapter. Hold the needle assembly between your thumb and forefinger, slide out the assembly towards the rear of the needle carrier assembly.



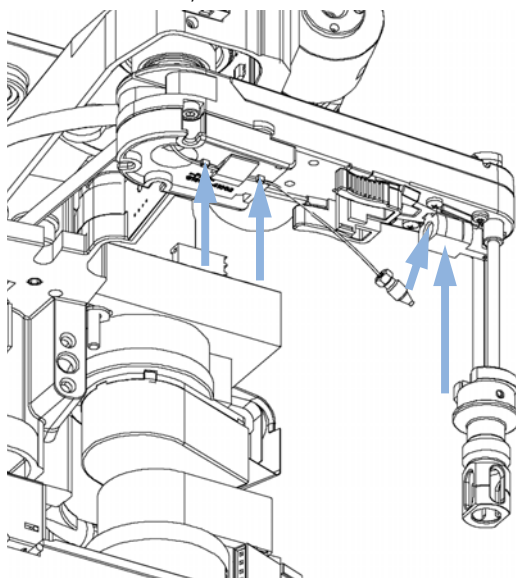
9 Maintenance

Exchanging the Needle Assembly

- 3** Insert the new needle assembly into the holder of the needle carrier assembly. Make sure to push it all the way to the front. Reassemble the wellplate adapter.



- 4** Use the 4 mm wrench and the 5/16" wrench for counter-holding to connect the valve to needle tubing to the needle. Reassemble the wellplate adapter. (Viewed from the bottom).



Next Steps:

- 5** **IMPORTANT:** After fixing the needle in the needle carrier clip the tubing into the guide on the bottom of the needle carrier assembly in case it slipped out of this guide. It is absolutely vital that the tubing is installed as described, to maximize the lifetime of the tubing.
- 6** Re-install the tray(s) in the tray base.
- 7** Start the instrument.
- 8** Close the front cover.

Exchanging the Needle/Capillary Carrier Assembly

When	If defective	
Tools required	p/n	Description
	8710-2438	Hex key 2.0 mm
Parts required	p/n	Description
	G1364-60022	Needle carrier assembly, analytical scale
Preparations	- Position the transport unit of the fraction collector in the Home position (see “Test Functions” on page 117). - Remove all installed trays from the tray base. - Position the transport unit of the fraction collector in the Change Parts position (see “Test Functions” on page 117) and turn OFF the instrument. - It might be more convenient to remove the needle from its carrier before unscrewing the needle tubing.	

WARNING

Personal injury

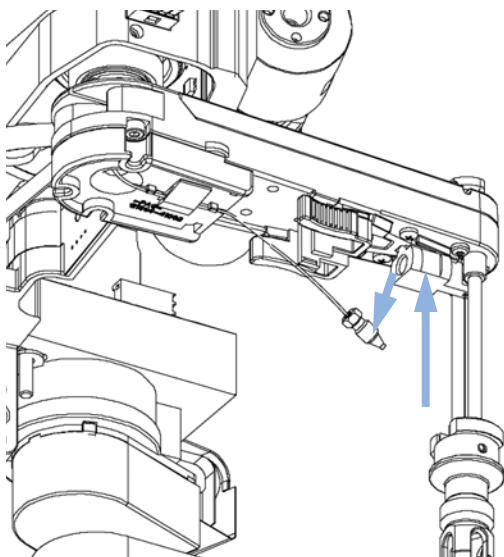
There is a risk of personal injury because the needle arm can move during fraction collector operation.

→ To avoid personal injury, keep fingers away from the needle area.

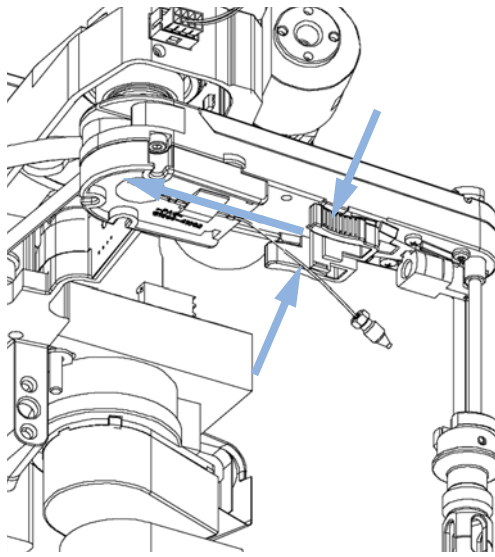
9 Maintenance

Exchanging the Needle/Capillary Carrier Assembly

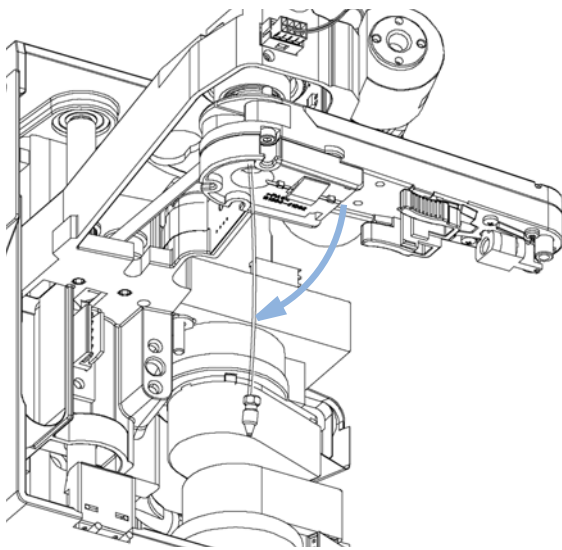
- 1** Use the 4 mm wrench and the 5/16" wrench for counter-holding to unscrew the valve-to-needle tubing from the needle assembly. (Viewed from the bottom).



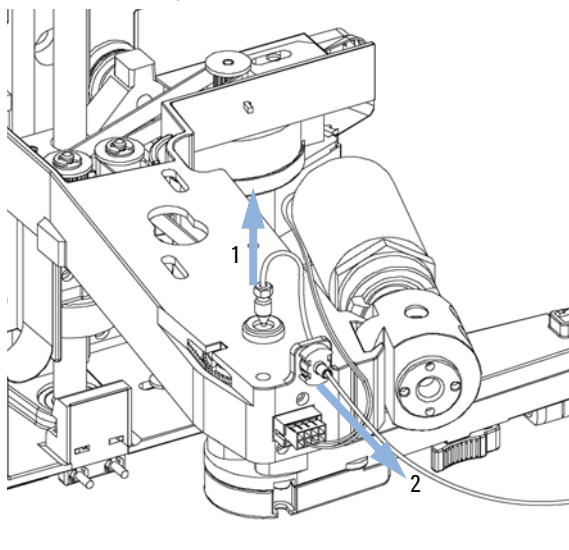
- 2** Hold the needle/capillary guiding assembly between your thumb and forefinger and slide out the assembly towards the rear of the needle carrier assembly.



- 3** Un-clip the tubing or capillary from the bottom of the needle carrier assembly.

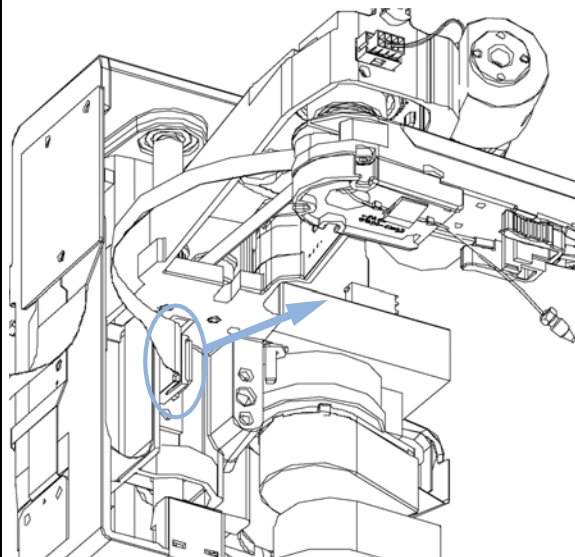


- 4** Slide the tubing or capillary through the hole in the carrier assembly (from bottom to top) and out of the holder in the z-arm assembly.

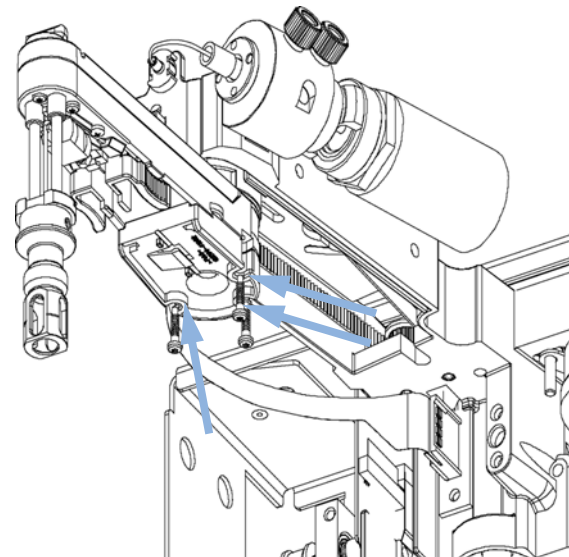


Exchanging the Needle/Capillary Carrier Assembly

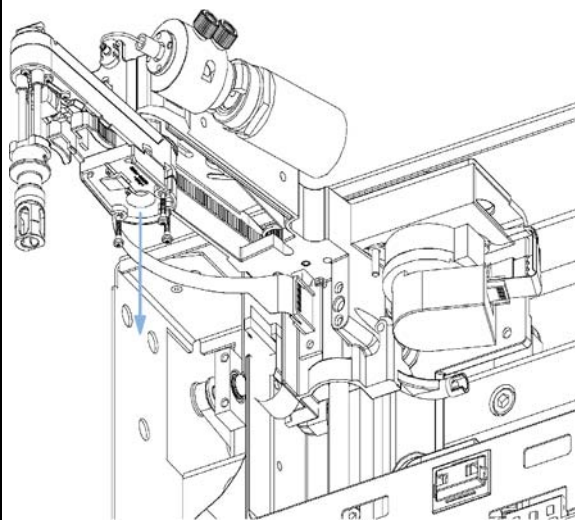
- 5** Unplug the blue connector of the brown flat band cable of the needle carrier assembly from its plug at the z-arm.



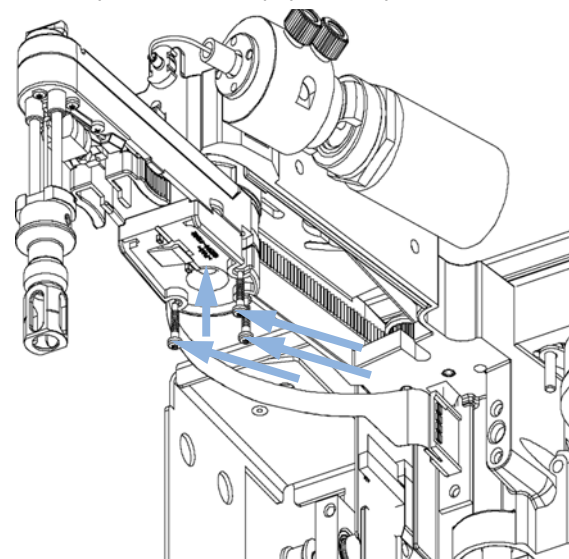
- 6** Locate the three holding screws of the needle carrier assembly and unscrew them (viewed from the bottom).



- 7** Slide out the (old) needle carrier assembly to the bottom.



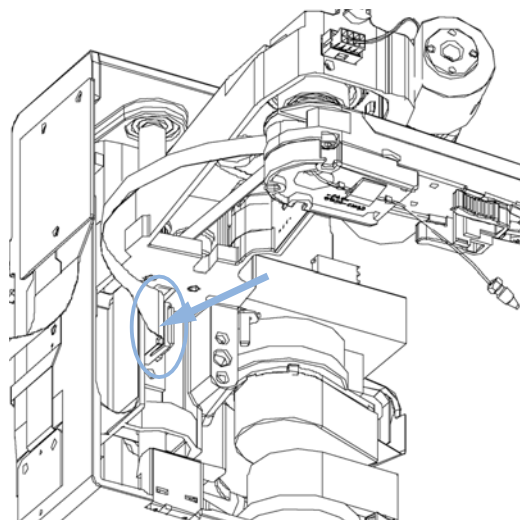
- 8** Install the new carrier assembly and secure it with its three holding screws. The carrier is held in position with a pin / slot - fit. It is only oriented and installed correctly, when pushed all the way up to the top.



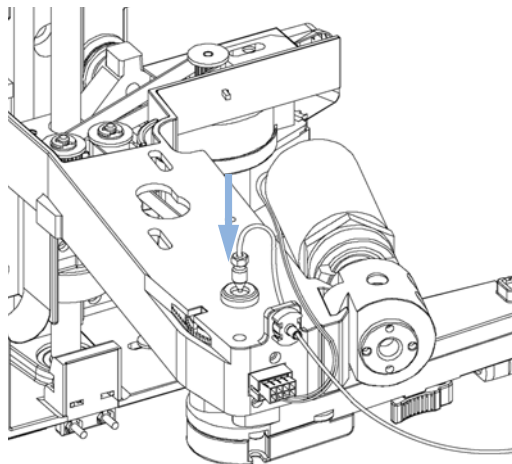
9 Maintenance

Exchanging the Needle/Capillary Carrier Assembly

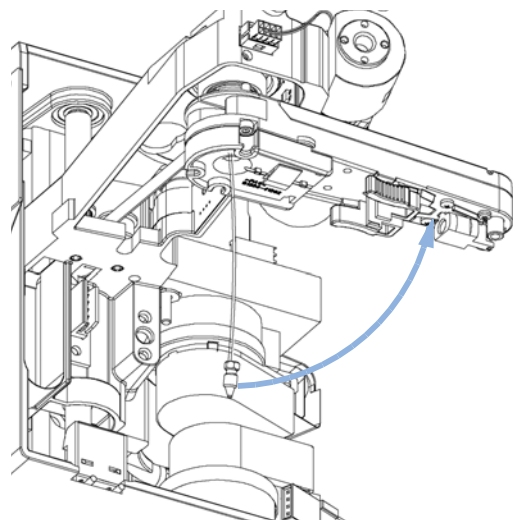
- 9** Plug in the blue connector of the brown flat band cable of the needle carrier assembly to its plug at the z-arm.



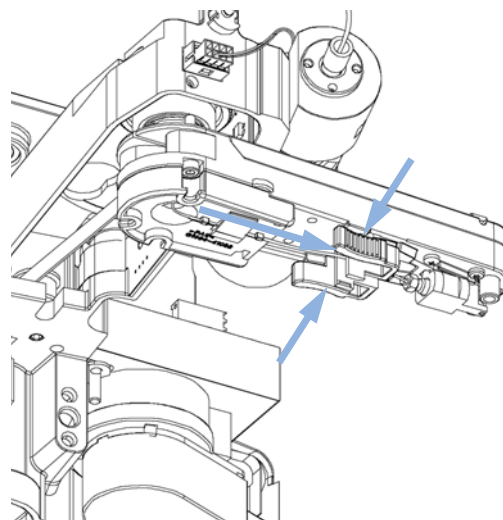
- 10** Slide the valve to needle tubing assembly through the hole in the z-arm and out on the bottom of the needle carrier assembly (top to bottom).



- 11** With the needle held underneath the needle carrier assembly and using the 4 mm wrench and the 5/16" wrench for counter-holding connect the valve to needle tubing to the needle. (Viewed from the bottom).

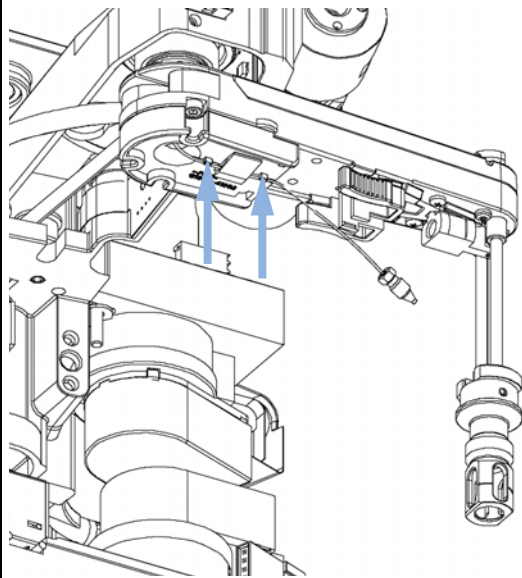


- 12** Insert the needle or capillary guiding assembly into the holder of the new needle carrier assembly. Make sure to push it all the way to the front (clicks into position).



Exchanging the Needle/Capillary Carrier Assembly

13 IMPORTANT: Clip the tubing or capillary into the guide on the bottom of the needle carrier assembly. It is absolutely vital that the tubing is installed as described, to maximize the lifetime of the tubing.

**Next Steps:**

14 IMPORTANT: After fixing the needle in the needle carrier clip the tubing into the guide on the bottom of the needle carrier assembly in case it slipped out of this guide. It is absolutely vital that the tubing is installed as described, to maximize the lifetime of the tubing.

15 Re-install the tray(s) in the tray base.

16 Start the instrument.

17 Close the front cover.

Exchanging the Diverter Valve

When If leaky or defective

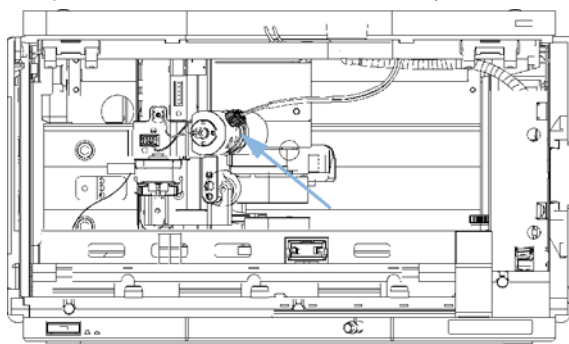
Tools required	p/n	Description
	8710-2438	Hex key 2.0 mm

Parts required	p/n	Description
	G5664-60901	3/2 Way valve assembly, bio-inert

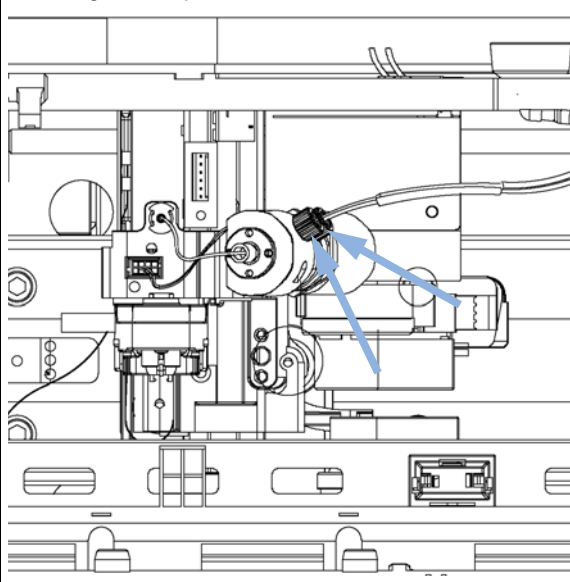
Preparations

- Position the transport unit of the fraction collector in the **Home** position (see “[Maintenance Functions](#)” on page 118).
- Remove all installed trays from the tray base.
- Turn OFF the instrument.

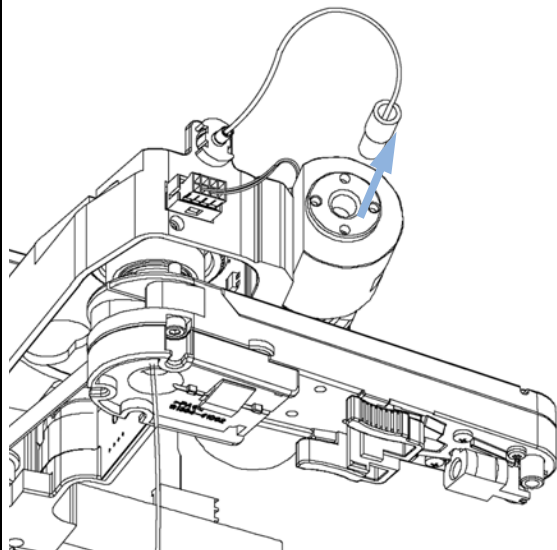
1 Locate the diverter valve with the finger tight fittings of the inlet / waste tubing assembly (the figure shows the open fraction collector seen from the front).



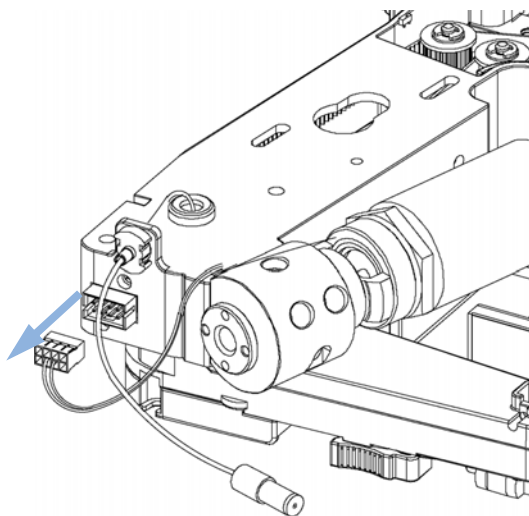
2 Unscrew the two finger-tight fittings of the inlet / waste tubing assembly at the diverter valve.



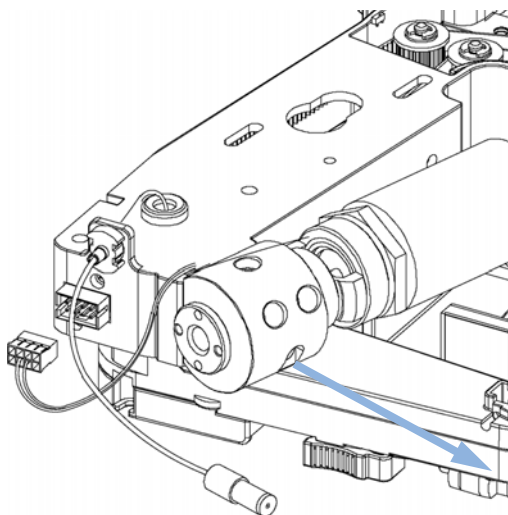
- 3** Unscrew the finger-tight fitting of the valve to needle tubing assembly at the diverter valve.



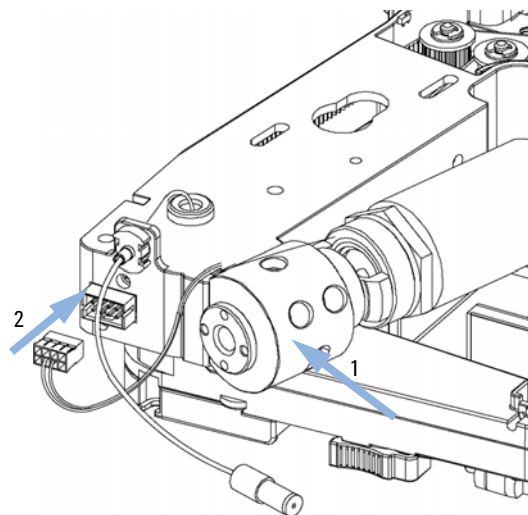
- 4** Disconnect the power cable of the diverter valve from the connector on the front of the z-arm.



- 5** Using the 2 mm hex key loosen the holding screw of the diverter valve and remove the diverter valve from the z-arm.



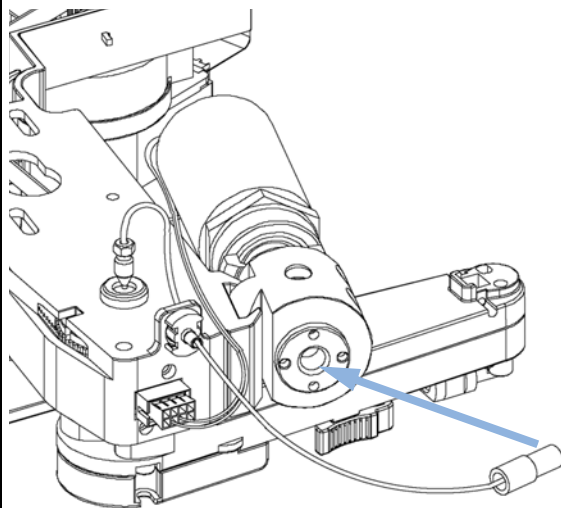
- 6** Install the new diverter valve to the z-arm with its holding screw. Plug in the new diverter valve power cable to the z-arm.



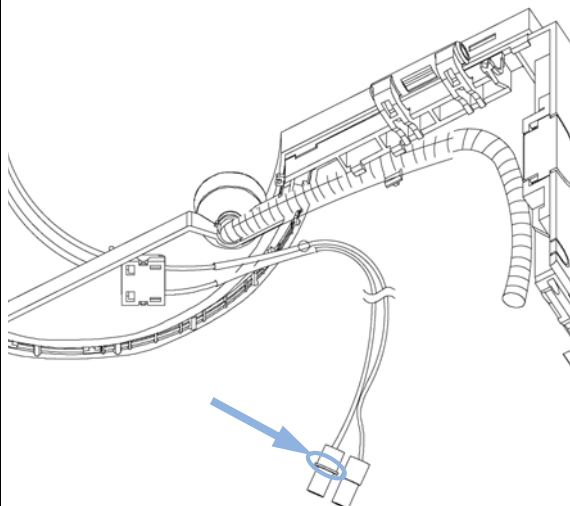
9 Maintenance

Exchanging the Diverter Valve

- 7** Install the valve to needle tubing at the new diverter valve.



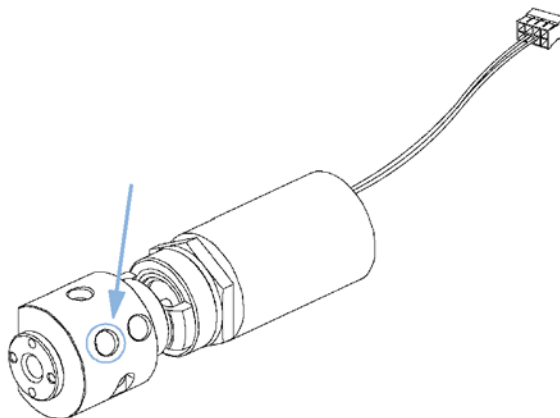
- 8** Connect the finger-tight fittings of the inlet / waste tubing assembly to the ports of the diverter valve. (View from the bottom)



NOTE

The tubings must not be bent up- or downwards. The cables must not be twisted.

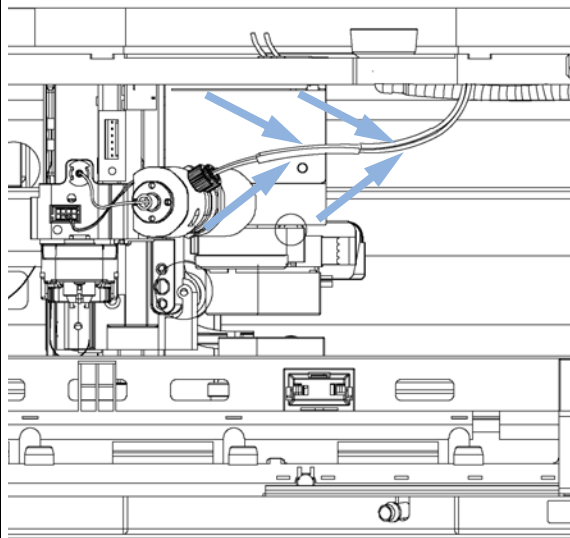
- 9** A color coded ring on one of the tubings and the valve body indicates, which cable belongs to which port.



NOTE

It is absolutely vital to connect these tubings as described, in order to maximize their lifetime and operating security.

10 IMPORTANT: The cables must run into the ports of the diverter valve in lines parallel to the horizon.



Next Steps:

11 Re-install the tray(s) in the tray base.

12 Start the instrument.

13 Close the front cover.

Exchanging the Internal Tray

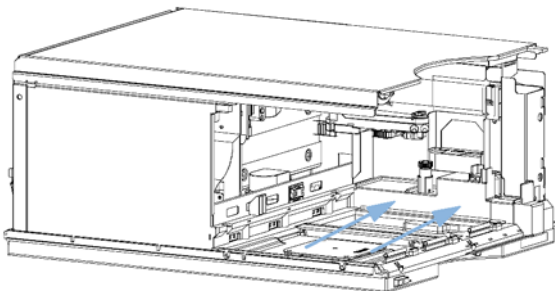
When If defective

Parts required	p/n	Description
	G1364-63114	Internal tray analytical scale

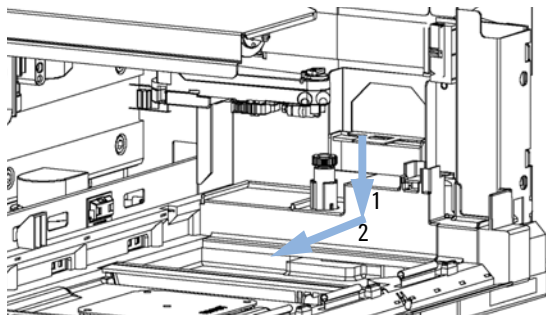
Preparations

- Position the transport unit of the fraction collector in the **Home** position (see [“Maintenance Functions”](#) on page 118).
- Remove all installed trays from the tray base.
- Turn OFF the instrument.

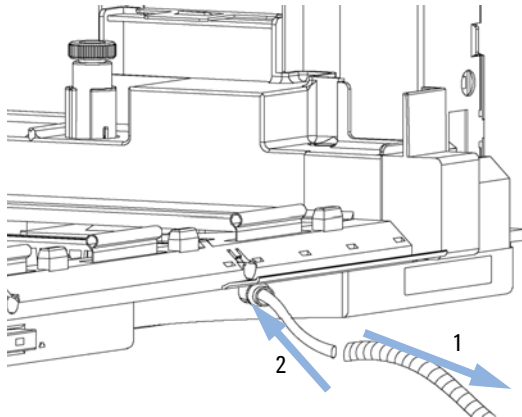
1 Locate the internal tray assembly with the rinse funnel and flow delay sensor in the bottom of the right front corner of the instrument.



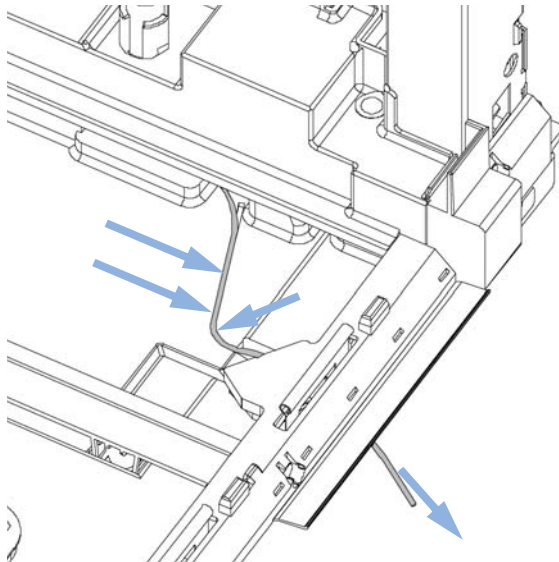
2 Remove the internal tray by pushing down the plastic holder that holds it in position underneath the metal leash (1) and sliding the tray to the left at the same time (2).



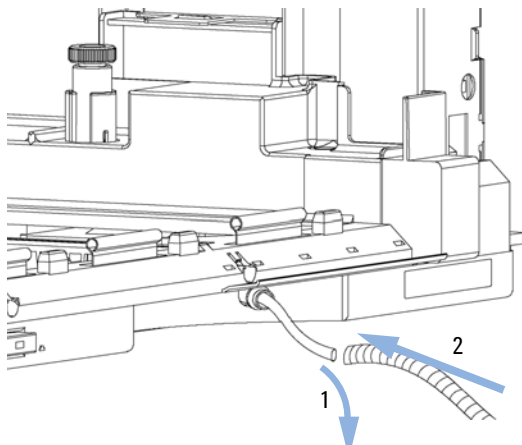
- 3** Remove the corrugated waste tubing from the front of the instrument (1) and slide the internal tray's waste tubing to the rear of the outlet (2) the before sliding the tray out.



- 4** Install the new tray by sliding it into position underneath the metal leash that holds it. The waste tubing from the internal tray should be guided as shown.



- 5** Make sure that the waste tubing is slid all the way through the outlet (1). Its end should be over the edge and below the level of the laboratory desk that the system stands on to avoid any back flow of solvent. Re-install the corrugated waste tubing (2).



Next Steps:

- 6** Re-install the tray(s) in the tray base.
- 7** Start the instrument.
- 8** Close the front cover.

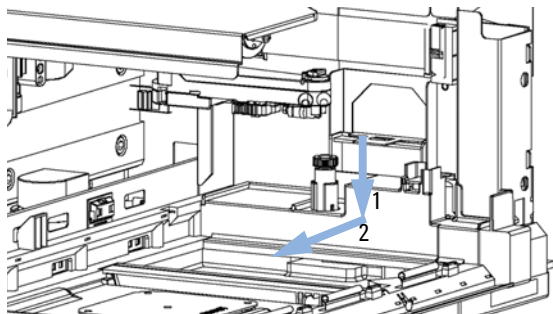
Repairing or Exchanging a Funnel of the Internal Tray

When When leaky or contaminated

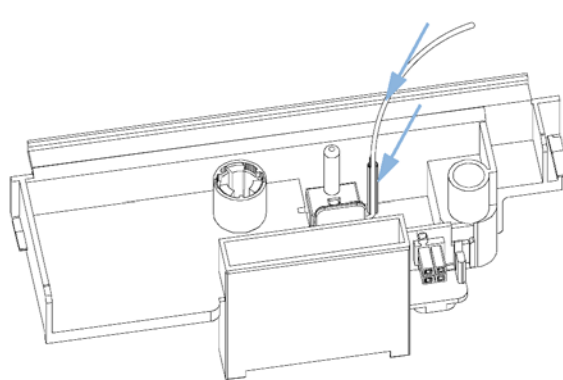
Parts required	p/n	Description
	G1364-68730	Funnel seal kit (pack of 10)
	G1364-43201	Funnel coupler
	G1364-86708	Waste tubing kit 0.5T (analytical scale)
	5022-2200	Funnel

- Preparations**
- Position the transport unit of the fraction collector in the **Home** position (see [“Maintenance Functions”](#) on page 118) and remove all installed trays from the tray base.
 - Turn OFF the instrument.

1 Remove the internal tray from the instrument, see [“Exchanging the Internal Tray”](#) on page 152.

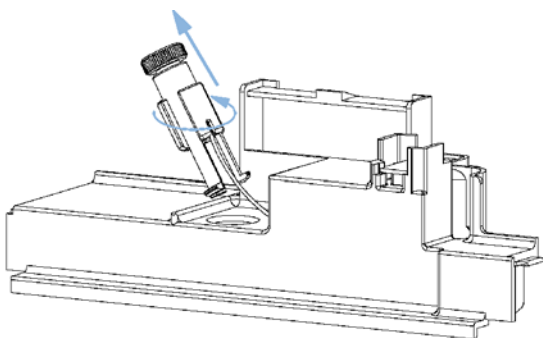


2 Turn the internal tray upside down and remove the funnel's waste tubing through the flow delay sensor.

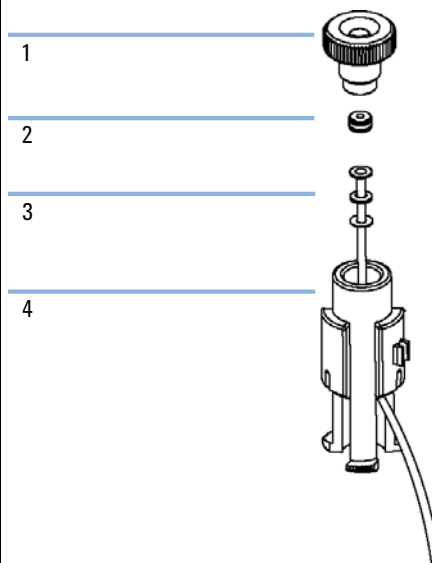


Repairing or Exchanging a Funnel of the Internal Tray

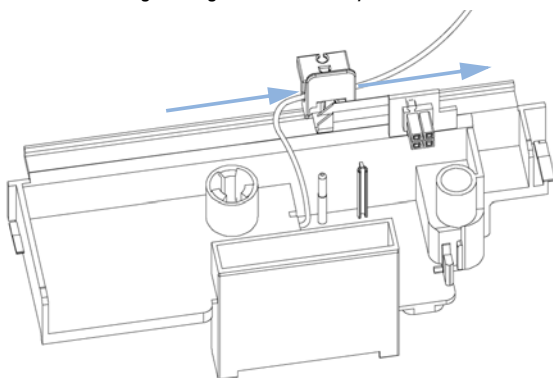
- 3** Remove the funnel assembly from the internal tray by screwing it counter clockwise and lifting it out.



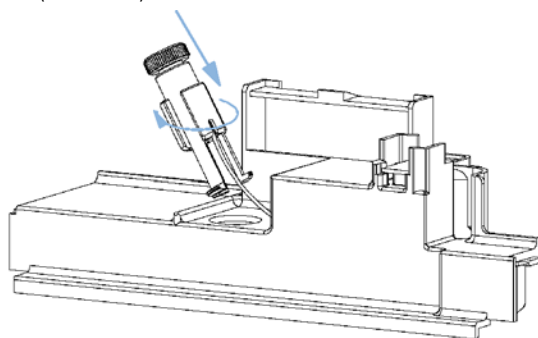
- 4** Exploded view for the parts of the internal tray. Once the funnel has been removed from the internal tray, it can be disassembled and defective parts can be replaced (funnel screw (1), seals (2), tubing (3) or funnel holder (4)).



- 5** Turn the internal tray upside down. Slide the funnel's waste tubing through the flow delay sensor.



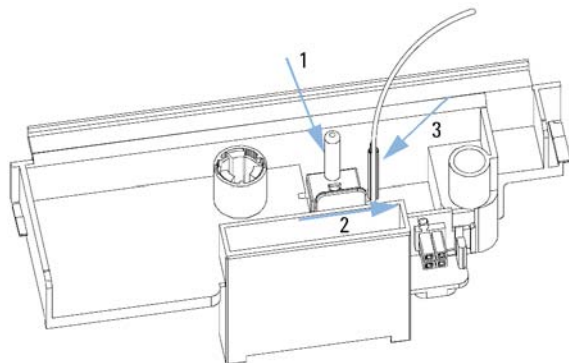
- 6** Insert the funnel into the tray and screw it tight (clockwise).



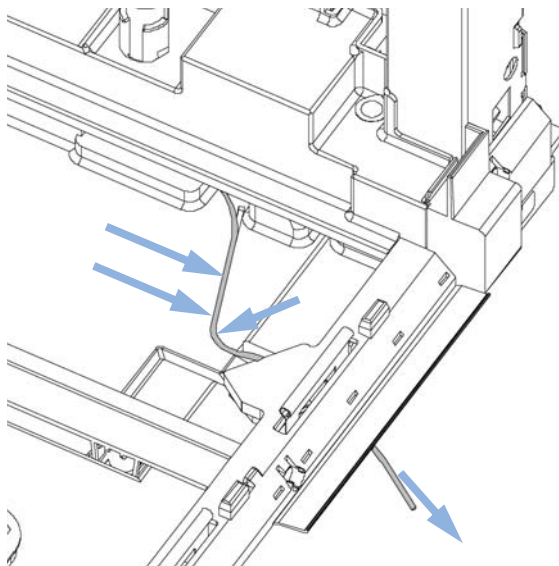
9 Maintenance

Repairing or Exchanging a Funnel of the Internal Tray

- 7** Re-install the plug that holds the flow delay sensor (1), pull the waste tubing tight through the flow delay sensor (2) and mount the waste tubing (3).



- 8** Re-install the internal tray and make sure to route the waste tubing out through the outlet of the instrument, correctly (see [“Exchanging the Internal Tray”](#) on page 152).



Next Steps:

- 9** Re-install the tray(s) in the tray base.
10 Start the instrument.
11 Close the front cover.

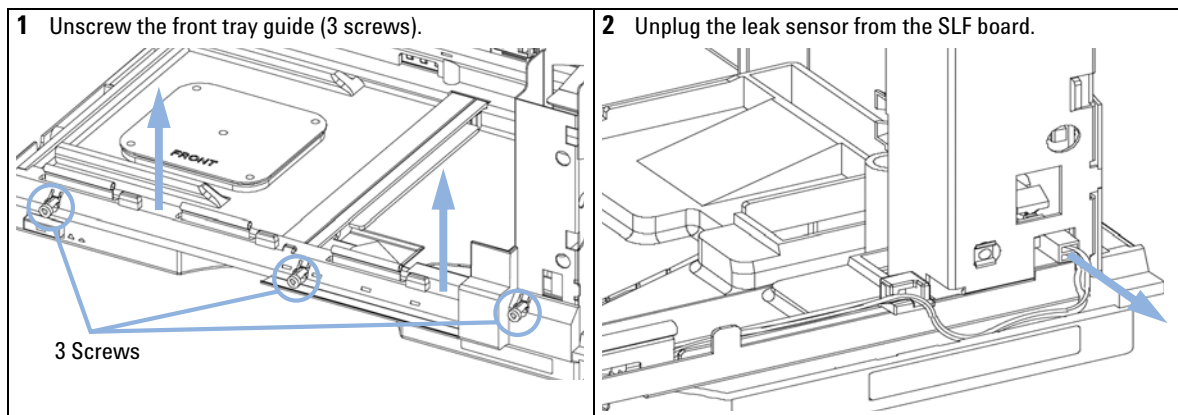
Exchanging the Leak Sensor

When If defective

Parts required	p/n	Description
	5061-3356	Leak sensor

Preparations

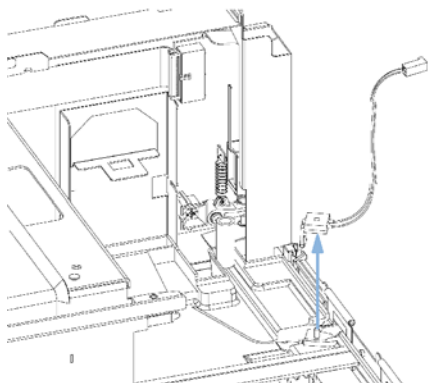
- Position the transport unit of the fraction collector in the **Home** position (see ["Maintenance Functions"](#) on page 118).
- Remove all installed trays from the tray base.
- Turn OFF the instrument.



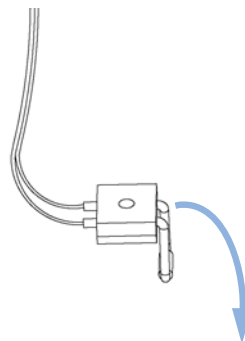
9 Maintenance

Exchanging the Leak Sensor

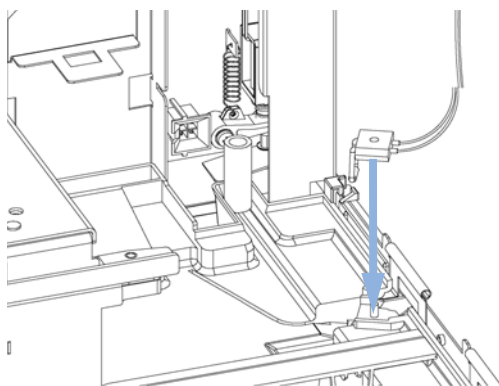
- 3** Notice the position of the leak sensor and pull out the old leak sensor and its cable from the instrument.



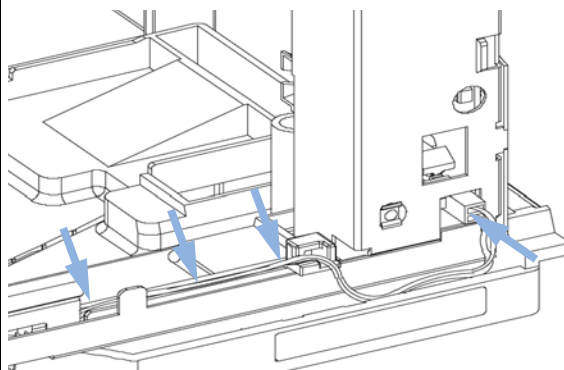
- 4** Bend the new leak sensor, so that it has a 90 ° angle to its plastic holder (as shown). The bend must be sharp enough that the sensor reaches all the way down to the bottom of the leak plane, but the cables must not be damaged!



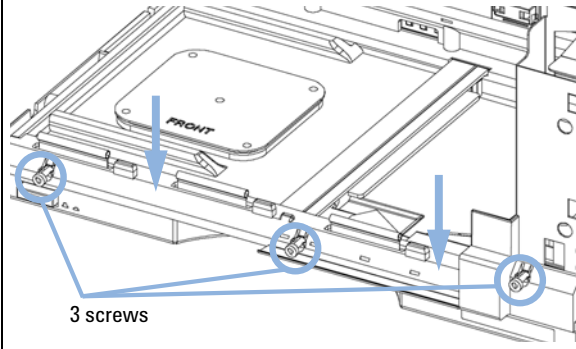
- 5** Install the new leak sensor into its position in the leak plane.



- 6** Route the leak sensor cable nicely into the recess at the front of the instrument. Plug the leak sensor into its connector on the SLF board.



- 7** Re-install the front tray guide (3 screws). Be careful not to pinch the new leak sensor's cable.



Next Steps:

- 8** Re-install the tray(s) in the tray base.
9 Start the instrument.
10 Close the front cover.

Exchanging or Installing the (Optional) Interface Board

When At installation, when defective or for all repairs inside the fraction collector.

Tools required **Description**
Screwdriver, flat-head

CAUTION

Electronic boards and components are sensitive to electrostatic discharge (ESD). ESD can damage electronic boards and components.

→ Be sure to hold the board by the edges, and do not touch the electrical components. Always use ESD protection (for example, an ESD wrist strap) when handling electronic boards and components.

-
- 1 Switch OFF the fraction collector at the main power switch.
 - 2 Disconnect cables from the interface board connectors.
 - 3 Loosen the screws. Slide out the interface board from the fraction collector.
 - 4 Install the interface board. Secure the screws.

Exchanging or Installing the (Optional) Interface Board

- 5 Reconnect the cables to the board connectors.

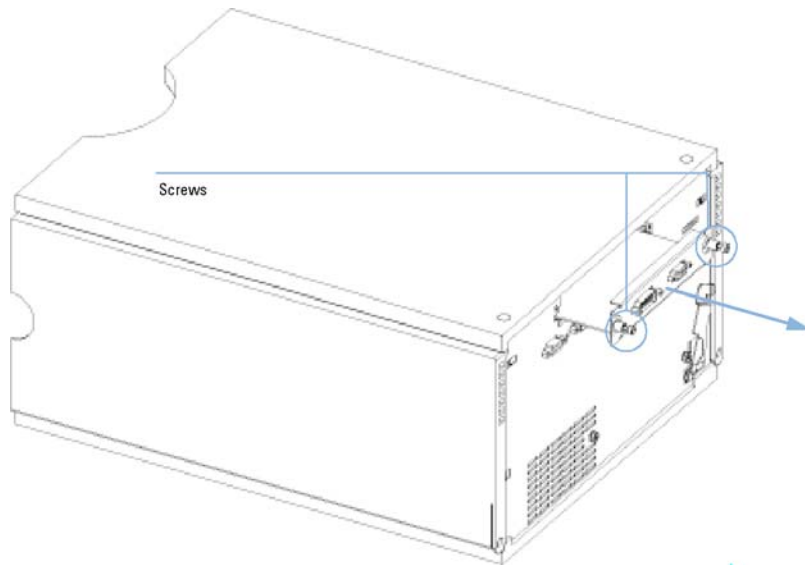


Figure 29 Exchanging the Interface Board

Replacing Module Firmware

When

The installation of newer firmware might be necessary

- if a newer version solves problems of older versions or
- to keep all systems on the same (validated) revision.

The installation of older firmware might be necessary

- to keep all systems on the same (validated) revision or
- if a new module with newer firmware is added to a system or
- if third part control software requires a special version.

Tools required

Description

LAN/RS-232 Firmware Update Tool

OR

Agilent Diagnostic Software

OR

Instant Pilot G4208A

(only if supported by module)

Parts required

Description

- 1 Firmware, tools and documentation from Agilent web site

Preparations

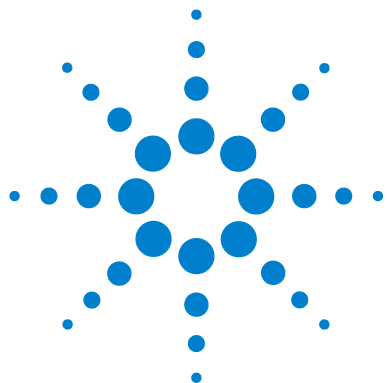
Read update documentation provided with the Firmware Update Tool.

To upgrade/downgrade the module's firmware carry out the following steps:

- 1 Download the required module firmware, the latest LAN/RS-232 FW Update Tool and the documentation from the Agilent web.
 - http://www.chem.agilent.com/scripts/cag_firmware.asp.
- 2 For loading the firmware into the module follow the instructions in the documentation.

Module Specific Information

There is no specific information for this module.



10 Parts for Maintenance

Fraction Collector Main Assemblies	164
Supported Trays for a Fraction Collector	166
List of Recommended Vials and Caps	167
Recommended Plates and Closing Mats	170
Fraction Collector Accessory Kit	172
Fraction Collector Thermostat	173
Needle Assembly	174
Diverter-Valve Assembly	175
Tubing Kits	176
Internal Tray Assembly	177

This chapter provides information on parts for maintenance.



Fraction Collector Main Assemblies

Item	p/n	Description
1	G5664-60901	3/2 Way valve assembly, bio-inert
2	G1364-63114	Internal tray analytical scale
3	G1367-47200	Plug channel
4	G1329-43200	Adapter air channel
5	G5667-87200	Needle assembly (bio-inert)
6	G1364-60022	Needle carrier assembly, analytical scale
7	G1351-68701	Interface board (BCD) with external contacts and BCD outputs (OPTIONAL)



For bio-inert modules use bio-inert parts only!

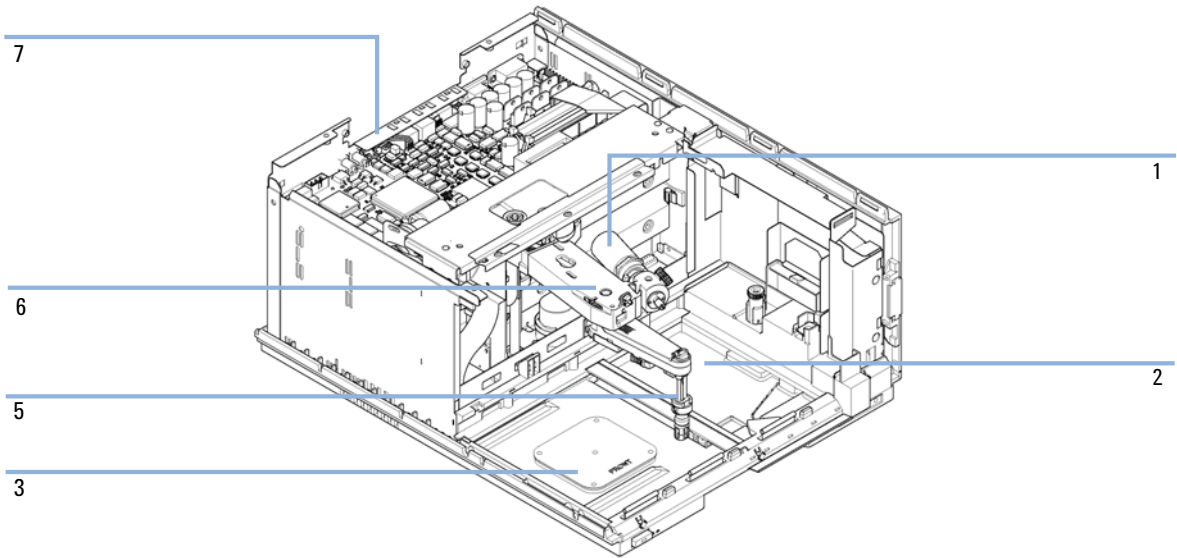


Figure 30 Fraction Collector Main Assemblies

Supported Trays for a Fraction Collector

For more information on trays, well-plates and collecting tubes for an LC Fraction Collector, refer to the Agilent Web site:

- <http://www.chem.agilent.com/en-US/Products/columns-supplies/instrument-parts/lc/pages/gp10885.aspx>
- <http://www.chem.agilent.com/en-US/Products/columns-supplies/generalchromatography/pages/gp10886.aspx>

List of Recommended Vials and Caps

Crimp Top Vials

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5181-3375	Crimp Top Vial, 2 mL, clear glass, 100/Pack
5183-4491	Crimp Top Vial, 2 mL, clear glass, 1000/Pack
5182-0543	Crimp Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-4492	Crimp Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-4494	Crimp Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)
5182-3376	Crimp Top Vial, 2 mL, amber glass, write-on spot, 100/Pack
5183-4493	Crimp Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-4495	Crimp Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)

SnapTop Vials

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5182-0544	Snap Top Vial, 2 mL, clear glass, 100/Pack
5183-4504	Snap Top Vial, 2 mL, clear glass, 1000/Pack
5183-4507	Snap Top Vial, 2 mL, clear glass, 100/Pack (silanized)
5182-0546	Snap Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-4505	Snap Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-4508	Snap Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)

10 Parts for Maintenance

List of Recommended Vials and Caps

p/n	Description
5182-0545	Snap Top Vial, 2 mL, amber glass, write-on spot, 100/Pack
5183-4506	Snap Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-4509	Snap Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)

Screw Top Vials

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5182-0714	Screw Cap Vials, 2 mL, clear glass, 100/Pack
5183-2067	Screw Top Vial, 2 mL, clear glass, 1000/Pack
5183-2070	Screw Top Vial, 2 mL, clear glass, 100/Pack (silanized)
5182-0715	Screw Top Vial, 2 mL, clear glass, write-on spot, 100/Pack
5183-2068	Screw Top Vial, 2 mL, clear glass, write-on spot, 1000/Pack
5183-2071	Screw Top Vial, 2 mL, clear glass, write-on spot, 100/Pack (silanized)
5182-0716	Screw Cap Vial, 2 mL, amber glass, write-on spot, 100/pk
5183-2069	Screw Top Vial, 2 mL, amber glass, write-on spot, 1000/Pack
5183-2072	Screw Top Vial, 2 mL, amber glass, write-on spot, 100/Pack (silanized)

Crime Caps

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5181-1210	Crimp Cap, silver aluminum, septum (clear PTFE/red rubber), 100/Pack
5183-4498	Crimp Cap, silver aluminum, septum (clear PTFE/red rubber), 1000/Pack
5181-1215	Crimp Cap, blue aluminum, septum (clear PTFE/red rubber), 100/Pack

p/n	Description
5181-1216	Crimp Cap, green aluminum, septum (clear PTFE/red rubber), 100/Pack
5181-1217	Crimp Cap, red aluminum, septum (clear PTFE/red rubber), 100/Pack

Snap Caps

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5182-0550	Snap Cap, clear polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3458	Snap Cap, blue polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3457	Snap Cap, green polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-3459	Snap Cap, red polypropylene, septum (clear PTFE/red rubber), 100/Pack

Screw Caps

NOTE

Caps for Use with the Analytical Scale Fraction Collector, only!

p/n	Description
5182-0717	Screw Cap, blue polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0718	Screw Cap, green polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0719	Screw Cap, red polypropylene, septum (clear PTFE/red rubber), 100/Pack
5182-0720	Screw Cap, blue polypropylene, septum (clear PTFE/silicone), 100/Pack
5182-0721	Screw Cap, green polypropylene, septum (clear PTFE/silicone), 100/Pack
5182-0722	Screw Cap, red polypropylene, septum (clear PTFE/silicone), 100/Pack

Recommended Plates and Closing Mats

Warnings

WARNING

Explosive gas mixtures

There is a risk of explosive gas mixtures in the instrument if flammable solvents are used.

- Cover the plates.
 - Remove the plates from the fraction collector after turning it OFF.
 - Only use solvents with a flash point higher than 200 °C.
-

WARNING

Contamination with adhesives

Closing mats with adhesives can give some contamination in the system. The adhesive is soluble in most of the solvents used in HPLC.

- In general do not use closing mats with adhesive. The fraction collector has no prepunch needle, therefore the adhesive will clog the needle after several injections.
-

List of Recommended Plates and Closing Mats

Recommended Plates and Closing Mats (Std. Well Plates and Closing Mats for Use with the Analytical Scale Fraction Collector, only!)

p/n	Description
5042-1386	96 well plate 0.5 ml, PP (pack of 10)
5042-1385	96 well plate 0.5 ml, PP (pack of 120)
5042-6454	96DeepAgilent31mm
5065-4402	96CappedAgilent
5188-5321	Glass inserts, 0.35 ml, 1000/Pack
5042-1388	384Agilent
5042-8502	96Agilent conical
G2255-68700	Vial plate for 54 x 2 mL vials (6/pk)
5022-6539	Vial plate for 15 x 6 mL vials (1/pk)
5022-6538	Vial plate for 27 Eppendorf tubes (1/pk)
5042-1389	Closing mat for all 96 Agilent plates

NOTE

Only one type of well-plates can be used at a time in one tray.

Fraction Collector Accessory Kit

p/n	Description
G1364-68755	Accessory Kit

p/n	Description
0100-1711	1/16" x 10-32 Male Nut, PEEK
0100-1856	3/16 ID Tube Barb Y Ftg PP/OBS
5042-6421	Micro Clamp
5042-6472	Clamp Waste Tube
5063-6527	Tubing assembly, i.d. 6 mm, o.d. 9 mm, 1.2 m (to waste)
5181-1519	CAN cable, Agilent module to module, 1 m
8710-2476	HEX-ALLEN-KEY 2.0-MM-A/F
G1313-44512	Halftray for 40 x 2 mL vials
G1313-44513	Halftray for 15 x 6 mL vials
G1329-43200	Adapter air channel

Fraction Collector Thermostat

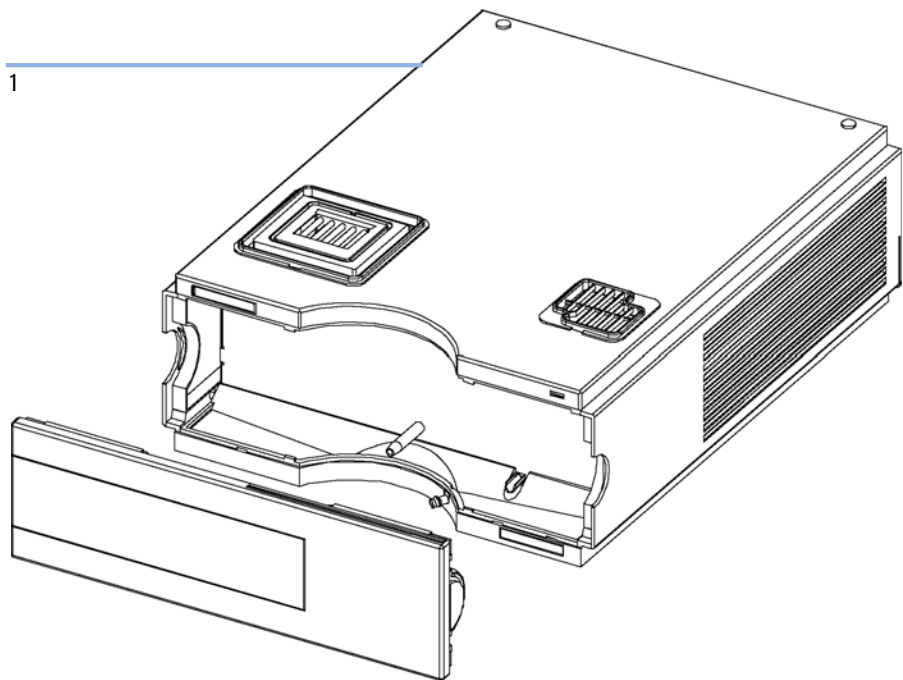


Figure 31 Fraction Collector Thermostat

Item	p/n	Description
1	G1330-69030	FRC thermostat, exchange assembly

Needle Assembly

p/n	Description
G5667-87200	Needle assembly (bio-inert)



For bio-inert modules use bio-inert parts only!

Diverter-Valve Assembly

Item	p/n	Description
1	G5664-60901	3/2 Way valve assembly, bio-inert
2	0515-1211	PIN screw



For bio-inert modules use bio-inert parts only!

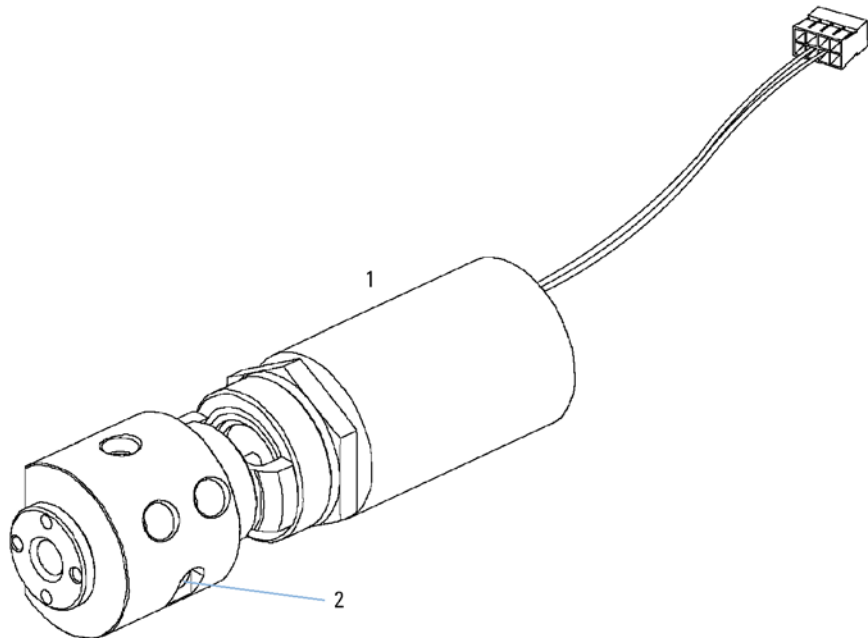


Figure 32 Diverter-Valve Assembly

Tubing Kits

Item	p/n	Description
	G5664-68712	Analytical tubing kit 0.25 mm i.d. PTFE-ESD
1	G5664-86706	Inlet/waste tubing assembly analytical scale 0.25 mm i.d.
2	G5664-86703	Diverter Valve to needle tubing analytical scale 0.25 mm i.d.
3	0100-1516	Fitting male PEEK, 2/pk

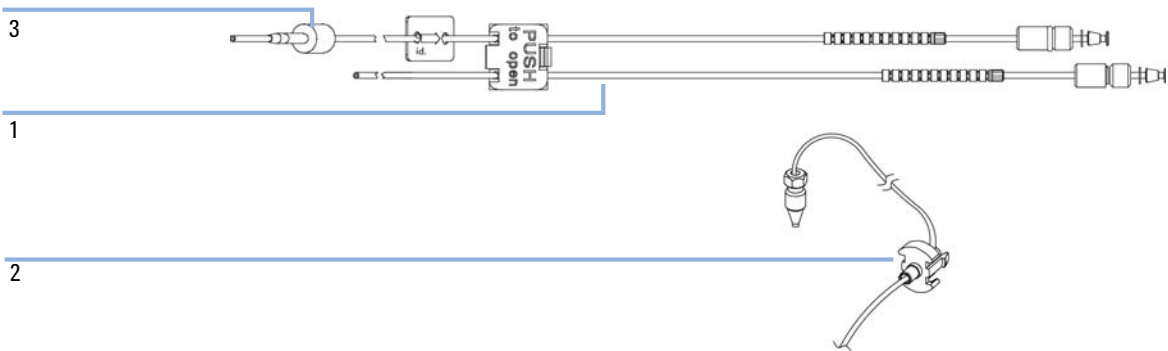


Figure 33 Tubing kits

Internal Tray Assembly

Item	p/n	Description
1	G1364-63114	Internal tray analytical scale
2	5022-2200	Funnel
3	G1364-68730	Funnel seal kit (pack of 10)
4	G1364-86708	Waste tubing kit 0.5T (analytical scale)
5	G1364-43201	Funnel coupler
		Flow delay sensor (not shown)

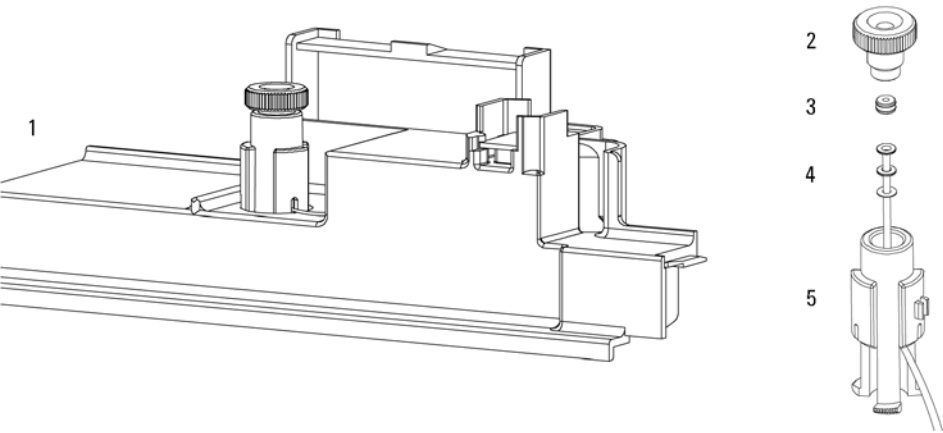


Figure 34 Internal Tray Assembly (Analytical Scale)

10 **Parts for Maintenance**
Internal Tray Assembly



11 Identifying Cables

Cable Overview	180
Analog Cables	182
Remote Cables	184
BCD Cables	187
CAN/LAN Cables	189
External Contact Cable	190
RS-232 Cable Kit	191

This chapter provides information on cables used with the Agilent modules.



Cable Overview

NOTE

Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.

Analog cables

p/n	Description
35900-60750	Agilent module to 3394/6 integrators
35900-60750	Agilent 35900A A/D converter
01046-60105	Analog cable (BNC to general purpose, spade lugs)

Remote cables

p/n	Description
03394-60600	Agilent module to 3396A Series I integrators 3396 Series II / 3395A integrator, see details in section “Remote Cables” on page 184
03396-61010	Agilent module to 3396 Series III / 3395B integrators
5061-3378	Remote Cable
01046-60201	Agilent module to general purpose

BCD cables

p/n	Description
03396-60560	Agilent module to 3396 integrators
G1351-81600	Agilent module to general purpose

CAN cables

p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

LAN cables

p/n	Description
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5023-0202	Twisted pair network cable, shielded, 7 m (for point to point connection)

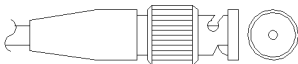
External Contact Cable

p/n	Description
G1103-61611	External contact cable - Agilent module interface board to general purposes

RS-232 cables

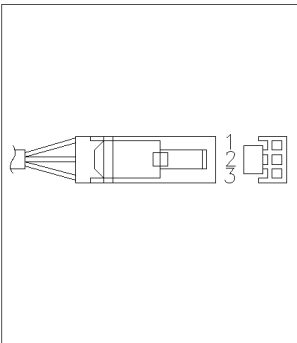
p/n	Description
G1530-60600	RS-232 cable, 2 m
RS232-61600	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It's also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 3-2, 4-6, 5-5, 6-4, 7-8, 8-7, 9-9.
5181-1561	RS-232 cable, 8 m

Analog Cables



One end of these cables provides a BNC connector to be connected to Agilent modules. The other end depends on the instrument to which connection is being made.

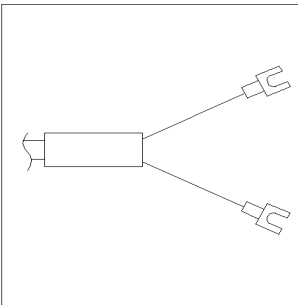
Agilent Module to 3394/6 Integrators

p/n 35900-60750	Pin 3394/6	Pin Agilent module	Signal Name
	1		Not connected
	2	Shield	Analog -
	3	Center	Analog +

Agilent Module to BNC Connector

p/n 8120-1840	Pin BNC	Pin Agilent module	Signal Name
	Shield	Shield	Analog -
	Center	Center	Analog +

Agilent Module to General Purpose

p/n 01046-60105	Pin	Pin Agilent module	Signal Name
	1		Not connected
	2	Black	Analog -
	3	Red	Analog +

Remote Cables



One end of these cables provides a Agilent Technologies APG (Analytical Products Group) remote connector to be connected to Agilent modules. The other end depends on the instrument to be connected to.

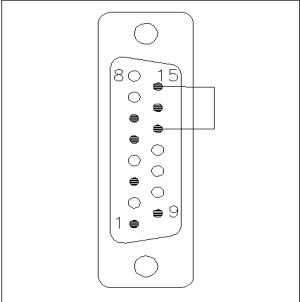
Agilent Module to 3396A Integrators

p/n 03394-60600	Pin 3396A	Pin Agilent module	Signal Name	Active (TTL)
	9	1 - White	Digital ground	
	NC	2 - Brown	Prepare run	Low
	3	3 - Gray	Start	Low
	NC	4 - Blue	Shut down	Low
	NC	5 - Pink	Not connected	
	NC	6 - Yellow	Power on	High
	5,14	7 - Red	Ready	High
	1	8 - Green	Stop	Low
	NC	9 - Black	Start request	Low
	13, 15		Not connected	

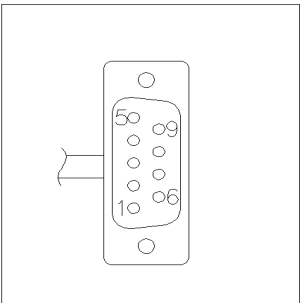
Agilent Module to 3396 Series II / 3395A Integrators

Use the cable Agilent module to 3396A Series I integrators (p/n 03394-60600) and cut pin #5 on the integrator side. Otherwise the integrator prints START; not ready.

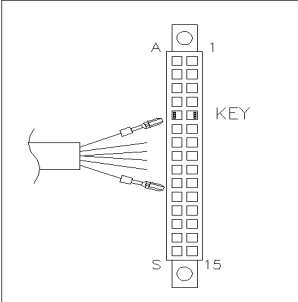
Agilent Module to 3396 Series III / 3395B Integrators

p/n 03396-61010	Pin 33XX	Pin Agilent module	Signal Name	Active (TTL)
	9	1 - White	Digital ground	
	NC	2 - Brown	Prepare run	Low
	3	3 - Gray	Start	Low
	NC	4 - Blue	Shut down	Low
	NC	5 - Pink	Not connected	
	NC	6 - Yellow	Power on	High
	14	7 - Red	Ready	High
	4	8 - Green	Stop	Low
	NC	9 - Black	Start request	Low
	13, 15		Not connected	

Agilent Module to Agilent 35900 A/D Converters

p/n 5061-3378	Pin 35900 A/D	Pin Agilent module	Signal Name	Active (TTL)
	1 - White	1 - White	Digital ground	
	2 - Brown	2 - Brown	Prepare run	Low
	3 - Gray	3 - Gray	Start	Low
	4 - Blue	4 - Blue	Shut down	Low
	5 - Pink	5 - Pink	Not connected	
	6 - Yellow	6 - Yellow	Power on	High
	7 - Red	7 - Red	Ready	High
	8 - Green	8 - Green	Stop	Low
	9 - Black	9 - Black	Start request	Low

Agilent Module to General Purpose

p/n 01046-60201	Wire Color	Pin Agilent module	Signal Name	Active (TTL)
	White	1	Digital ground	
	Brown	2	Prepare run	Low
	Gray	3	Start	Low
	Blue	4	Shut down	Low
	Pink	5	Not connected	
	Yellow	6	Power on	High
	Red	7	Ready	High
	Green	8	Stop	Low
	Black	9	Start request	Low

BCD Cables

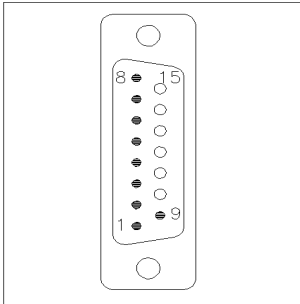


One end of these cables provides a 15-pin BCD connector to be connected to the Agilent modules. The other end depends on the instrument to be connected to

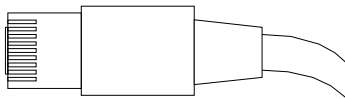
Agilent Module to General Purpose

p/n G1351-81600	Wire Color	Pin Agilent module	Signal Name	BCD Digit
	Green	1	BCD 5	20
	Violet	2	BCD 7	80
	Blue	3	BCD 6	40
	Yellow	4	BCD 4	10
	Black	5	BCD 0	1
	Orange	6	BCD 3	8
	Red	7	BCD 2	4
	Brown	8	BCD 1	2
	Gray	9	Digital ground	Gray
	Gray/pink	10	BCD 11	800
	Red/blue	11	BCD 10	400
	White/green	12	BCD 9	200
	Brown/green	13	BCD 8	100
	not connected	14		
	not connected	15	+ 5 V	Low

Agilent Module to 3396 Integrators

p/n 03396-60560	Pin 3396	Pin Agilent module	Signal Name	BCD Digit
	1	1	BCD 5	20
	2	2	BCD 7	80
	3	3	BCD 6	40
	4	4	BCD 4	10
	5	5	BCD0	1
	6	6	BCD 3	8
	7	7	BCD 2	4
	8	8	BCD 1	2
	9	9	Digital ground	
	NC	15	+ 5 V	Low

CAN/LAN Cables



Both ends of this cable provide a modular plug to be connected to Agilent modules CAN or LAN connectors.

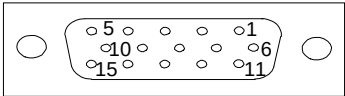
CAN Cables

p/n	Description
5181-1516	CAN cable, Agilent module to module, 0.5 m
5181-1519	CAN cable, Agilent module to module, 1 m

LAN Cables

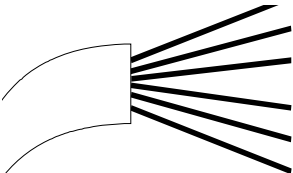
p/n	Description
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5023-0202	Twisted pair network cable, shielded, 7 m (for point to point connection)

External Contact Cable



One end of this cable provides a 15-pin plug to be connected to Agilent modules interface board. The other end is for general purpose.

Agilent Module Interface Board to general purposes

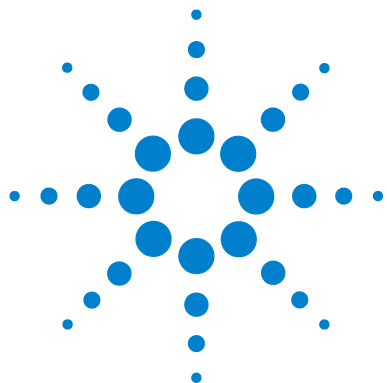
p/n G1103-61611	Color	Pin Agilent module	Signal Name
	White	1	EXT 1
	Brown	2	EXT 1
	Green	3	EXT 2
	Yellow	4	EXT 2
	Grey	5	EXT 3
	Pink	6	EXT 3
	Blue	7	EXT 4
	Red	8	EXT 4
	Black	9	Not connected
	Violet	10	Not connected
	Grey/pink	11	Not connected
	Red/blue	12	Not connected
	White/green	13	Not connected
	Brown/green	14	Not connected
	White/yellow	15	Not connected

RS-232 Cable Kit

p/n	Description
G1530-60600	RS-232 cable, 2 m
RS232-61600	RS-232 cable, 2.5 m Instrument to PC, 9-to-9 pin (female). This cable has special pin-out, and is not compatible with connecting printers and plotters. It's also called "Null Modem Cable" with full handshaking where the wiring is made between pins 1-1, 2-3, 3-2, 4-6, 5-5, 6-4, 7-8, 8-7, 9-9.
5181-1561	RS-232 cable, 8 m

11 Identifying Cables

RS-232 Cable Kit



12 Hardware Information

Electrical Connections	194
Rear view of the module	195
Interfaces	196
Overview Interfaces	198
Setting the 8-bit Configuration Switch (without On-board) LAN	202
Communication Settings for RS-232C	203
Special Settings	205
Optional Interface Boards	206

This chapter describes the module in more detail on hardware and electronics.



Electrical Connections

- The CAN bus is a serial bus with high speed data transfer. The two connectors for the CAN bus are used for internal module data transfer and synchronization.
- One analog output provides signals for integrators or data handling systems.
- The REMOTE connector may be used in combination with other analytical instruments from Agilent Technologies if you want to use features such as start, stop, common shut down, prepare, and so on.
- With the appropriate software, the RS-232C connector may be used to control the module from a computer through a RS-232C connection. This connector is activated and can be configured with the configuration switch.
- The power input socket accepts a line voltage of 100 – 240 VAC $\pm$ 10 % with a line frequency of 50 or 60 Hz. Maximum power consumption varies by module. There is no voltage selector on your module because the power supply has wide-ranging capability. There are no externally accessible fuses, because automatic electronic fuses are implemented in the power supply.

NOTE

Never use cables other than the ones supplied by Agilent Technologies to ensure proper functionality and compliance with safety or EMC regulations.

Rear view of the module

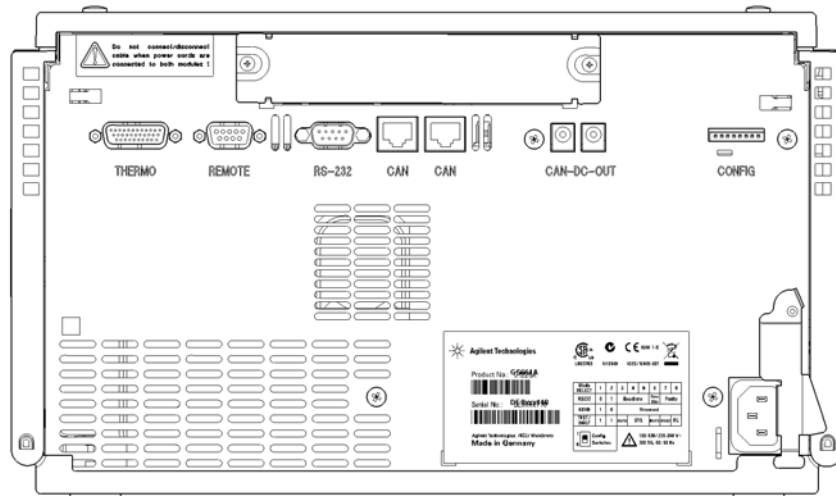


Figure 35 Fraction Collector Electrical Connections

Interfaces

The Agilent 1200 Infinity Series modules provide the following interfaces:

Table 17 Agilent 1200 Infinity Series Interfaces

Module	CAN	LAN/BCD (optional)	LAN (on-board)	RS-232	Analog	APG Remote	Special
Pumps							
G1310B Iso Pump	2	Yes	No	Yes	1	Yes	
G1311B Quat Pump							
G1311C Quat Pump VL							
G1312B Bin Pump							
G1312C Bin Pump VL							
1376A Cap Pump							
G2226A Nano Pump							
G5611A Bio-inert Quat Pump							
G4220A/B Bin Pump	2	No	Yes	Yes	No	Yes	
G1361A Prep Pump	2	Yes	No	Yes	No	Yes	CAN-DC- OUT for CAN slaves
Samplers							
G1329B ALS	2	Yes	No	Yes	No	Yes	THERMOSTAT for G1330B
G2260A Prep ALS							
G1364B FC-PS	2	Yes	No	Yes	No	Yes	THERMOSTAT for G1330B
G1364C FC-AS							CAN-DC- OUT for CAN slaves
G1364D FC- μ S							
G1367E HiP ALS							
G1377A HiP micro ALS							
G2258A DL ALS							
G5664A Bio-inert FC-AS							
G5667A Bio-inert Autosampler							
G4226A ALS	2	Yes	No	Yes	No	Yes	

Table 17 Agilent 1200 Infinity Series Interfaces

Module	CAN	LAN/BCD (optional)	LAN (on-board)	RS-232	Analog	APG Remote	Special
Detectors							
G1314B VWD VL G1314C VWD VL+	2	Yes	No	Yes	1	Yes	
G1314E/F VWD	2	No	Yes	Yes	1	Yes	
G4212A/B DAD	2	No	Yes	Yes	1	Yes	
G1315C DAD VL+ G1365C MWD G1315D DAD VL G1365D MWD VL	2	No	Yes	Yes	2	Yes	
G1321B FLD G1362A RID	2	Yes	No	Yes	1	Yes	
G4280A ELSD	No	No	No	Yes	Yes	Yes	EXT Contact AUTOZERO
Others							
G1170A Valve Drive	2	No	No	No	No	No	Requires a HOST module with on-board LAN (e.g. G4212A or G4220A with minimum firmware B.06.40 or C.06.40) or with additional G1369C LAN Card
G1316A/C TCC	2	No	No	Yes	No	Yes	
G1322A DEG	No	No	No	No	No	Yes	AUX
G1379B DEG	No	No	No	Yes	No	Yes	
G4225A DEG	No	No	No	Yes	No	Yes	
G4227A Flex Cube	2	No	No	No	No	No	
G4240A CHIP CUBE	2	Yes	No	Yes	No	Yes	CAN-DC- OUT for CAN slaves THERMOSTAT for G1330A/B (NOT USED)

NOTE

The detector (DAD/MWD/FLD/VWD/RID) is the preferred access point for control via LAN. The inter-module communication is done via CAN.

- CAN connectors as interface to other modules
- LAN connector as interface to the control software
- RS-232C as interface to a computer
- REMOTE connector as interface to other Agilent products
- Analog output connector(s) for signal output

Overview Interfaces

CAN

The CAN is inter-module communication interface. It is a 2-wire serial bus system supporting high speed data communication and real-time requirement.

LAN

The modules have either an interface slot for an LAN card (e.g. Agilent G1369B/C LAN Interface) or they have an on-board LAN interface (e.g. detectors G1315C/D DAD and G1365C/D MWD). This interface allows the control of the module/system via a PC with the appropriate control software.

NOTE

If an Agilent detector (DAD/MWD/FLD/VWD/RID) is in the system, the LAN should be connected to the DAD/MWD/FLD/VWD/RID (due to higher data load). If no Agilent detector is part of the system, the LAN interface should be installed in the pump or autosampler.

RS-232C (Serial)

The RS-232C connector is used to control the module from a computer through RS-232C connection, using the appropriate software. This connector can be configured with the configuration switch module at the rear of the module. Refer to *Communication Settings for RS-232C*.

NOTE

There is no configuration possible on main boards with on-board LAN. These are pre-configured for

- 19200 baud,
- 8 data bit with no parity and
- one start bit and one stop bit are always used (not selectable).

The RS-232C is designed as DCE (data communication equipment) with a 9-pin male SUB-D type connector. The pins are defined as:

Table 18 RS-232C Connection Table

Pin	Direction	Function
1	In	DCD
2	In	RxD
3	Out	TxD
4	Out	DTR
5		Ground
6	In	DSR
7	Out	RTS
8	In	CTS
9	In	RI

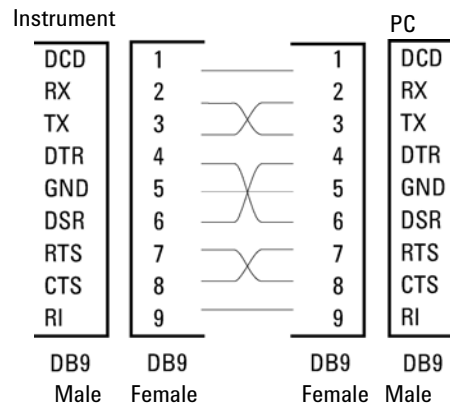


Figure 36 RS-232 Cable

Analog Signal Output

The analog signal output can be distributed to a recording device. For details refer to the description of the module's main board.

APG Remote

The APG Remote connector may be used in combination with other analytical instruments from Agilent Technologies if you want to use features as common shut down, prepare, and so on.

Remote control allows easy connection between single instruments or systems to ensure coordinated analysis with simple coupling requirements.

The subminiature D connector is used. The module provides one remote connector which is inputs/outputs (wired- or technique).

To provide maximum safety within a distributed analysis system, one line is dedicated to **SHUT DOWN** the system's critical parts in case any module detects a serious problem. To detect whether all participating modules are switched on or properly powered, one line is defined to summarize the **POWER ON** state of all connected modules. Control of analysis is maintained by signal readiness **READY** for next analysis, followed by **START** of run and optional **STOP** of run triggered on the respective lines. In addition **PREPARE** and **START REQUEST** may be issued. The signal levels are defined as:

- standard TTL levels (0 V is logic true, + 5.0 V is false),
- fan-out is 10,

- input load is 2.2 kOhm against + 5.0 V, and
- output are open collector type, inputs/outputs (wired- or technique).

NOTE

All common TTL circuits operate with a 5 V power supply. A TTL signal is defined as "low" or L when between 0 V and 0.8 V and "high" or H when between 2.0 V and 5.0 V (with respect to the ground terminal).

Table 19 Remote Signal Distribution

Pin	Signal	Description
1	DGND	Digital ground
2	PREPARE	(L) Request to prepare for analysis (for example, calibration, detector lamp on). Receiver is any module performing pre-analysis activities.
3	START	(L) Request to start run / timetable. Receiver is any module performing run-time controlled activities.
4	SHUT DOWN	(L) System has serious problem (for example, leak: stops pump). Receiver is any module capable to reduce safety risk.
5		Not used
6	POWER ON	(H) All modules connected to system are switched on. Receiver is any module relying on operation of others.
7	READY	(H) System is ready for next analysis. Receiver is any sequence controller.
8	STOP	(L) Request to reach system ready state as soon as possible (for example, stop run, abort or finish and stop injection). Receiver is any module performing run-time controlled activities.
9	START REQUEST	(L) Request to start injection cycle (for example, by start key on any module). Receiver is the autosampler.

Special Interfaces

Some modules have module specific interfaces/connectors. They are described in the module documentation.

Setting the 8-bit Configuration Switch (without On-board) LAN

The 8-bit configuration switch is located at the rear of the module.

This module does not have its own on-board LAN interface. It can be controlled through the LAN interface of another module, and a CAN connection to that module.

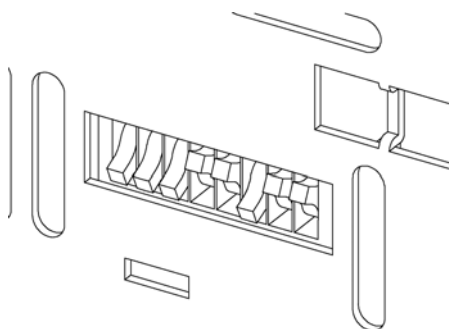


Figure 37 Configuration switch (settings depend on configured mode)

All modules without on-board LAN:

- default should be ALL DIPS DOWN (= best settings)
 - Bootp mode for LAN and
 - 19200 baud, 8 data bit / 1 stop bit with no parity for RS-232
- DIP 1 DOWN and DIP 2 UP allows special RS-232 settings
- for boot/test modes DIPS 1+2 must be UP plus required mode

NOTE

For normal operation use the default (best) settings.

Switch settings provide configuration parameters for serial communication protocol and instrument specific initialization procedures.

NOTE

With the introduction of the Agilent 1260 Infinity, all GPIB interfaces have been removed. The preferred communication is LAN.

NOTE

The following tables represent the configuration switch settings for the modules without on-board LAN only.

Table 20 8-bit Configuration Switch (without on-board LAN)

Mode Select	1	2	3	4	5	6	7	8
RS-232C	0	1	Baudrate			Data Bits	Parity	
Reserved	1	0	Reserved					
TEST/BOOT	1	1	RSVD	SYS		RSVD	RSVD	FC

NOTE

The LAN settings are done on the LAN Interface Card G1369B/C. Refer to the documentation provided with the card.

Communication Settings for RS-232C

The communication protocol used in the column compartment supports only hardware handshake (CTS/RTR).

Switches 1 in down and 2 in up position define that the RS-232C parameters will be changed. Once the change has been completed, the column instrument must be powered up again in order to store the values in the non-volatile memory.

Table 21 Communication Settings for RS-232C Communication (without on-board LAN)

Mode Select	1	2	3	4	5	6	7	8
RS-232C	0	1	Baudrate			Data Bits	Parity	

Use the following tables for selecting the setting which you want to use for RS-232C communication. The number 0 means that the switch is down and 1 means that the switch is up.

12 Hardware Information

Setting the 8-bit Configuration Switch (without On-board) LAN

Table 22 Baudrate Settings (without on-board LAN)

Switches			Baud Rate	Switches			Baud Rate
3	4	5		3	4	5	
0	0	0	9600	1	0	0	9600
0	0	1	1200	1	0	1	14400
0	1	0	2400	1	1	0	19200
0	1	1	4800	1	1	1	38400

Table 23 Data Bit Settings (without on-board LAN)

Switch 6	Data Word Size
0	7 Bit Communication
1	8 Bit Communication

Table 24 Parity Settings (without on-board LAN)

Switches		Parity
7	8	
0	0	No Parity
0	1	Odd Parity
1	1	Even Parity

One start bit and one stop bit are always used (not selectable).

Per default, the module will turn into 19200 baud, 8 data bit with no parity.

Special Settings

The special settings are required for specific actions (normally in a service case).

Boot-Resident

Firmware update procedures may require this mode in case of firmware loading errors (main firmware part).

If you use the following switch settings and power the instrument up again, the instrument firmware stays in the resident mode. It is not operable as a module. It only uses basic functions of the operating system for example, for communication. In this mode the main firmware can be loaded (using update utilities).

Table 25 Boot Resident Settings (without on-board LAN)

Mode Select	SW1	SW2	SW3	SW4	SW5	SW6	SW7	SW8
TEST/BOOT	1	1	0	0	1	0	0	0

Forced Cold Start

A forced cold start can be used to bring the module into a defined mode with default parameter settings.

CAUTION

Loss of data

Forced cold start erases all methods and data stored in the non-volatile memory. Exceptions are calibration settings, diagnosis and repair log books which will not be erased.

→ Save your methods and data before executing a forced cold start.

If you use the following switch settings and power the instrument up again, a forced cold start has been completed.

Table 26 Forced Cold Start Settings (without on-board LAN)

Mode Select	SW1	SW2	SW3	SW4	SW5	SW6	SW7	SW8
TEST/BOOT	1	1	0	0	1	0	0	1

Optional Interface Boards

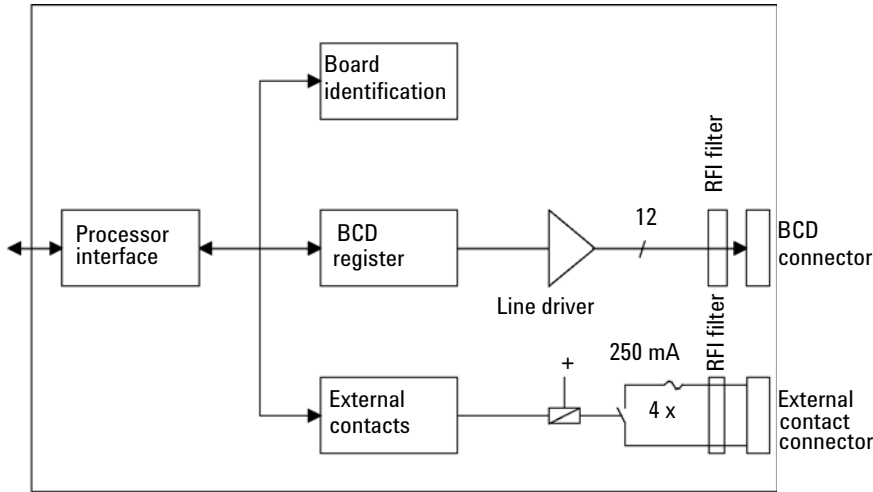
BCD / External Contact Board

The Agilent 1200 Infinity Series modules have one optional board slot that allows to add an interface board to the modules. Some modules do not have this interface slot. Refer to “[Interfaces](#)” on page 196 for details.

Optional Interface Boards

p/n	Description
G1351-68701	Interface board (BCD) with external contacts and BCD outputs
2110-0004	Fuse for BCD board, 250 mA

The BCD board provides a BCD output for the bottle number of the Agilent 1200 Series autosampler and four external contacts. The external contact closure contacts are relay contacts. The maximum settings are: 30 V (AC/DC); 250 mA (fused).



There are general purpose cables available to connect the BCD output, see “[BCD Cables](#)” on page 187 and the external outputs, see “[External Contact Cable](#)” on page 190 to external devices.

Table 27 Detailed connector layout (1200)

Pin	Signal name	BCD digit
1	BCD 5	20
2	BCD 7	80
3	BCD 6	40
4	BCD 4	10
5	BCD 0	1
6	BCD 3	8
7	BCD 2	4
8	BCD 1	2
9	Digital ground	
10	BCD 11	800
11	BCD 10	400
12	BCD 9	200
13	BCD 8	100
15	+5V	Low

LAN Communication Interface Board

The Agilent modules have one optional board slot that allows to add an interface board to the modules. Some modules do not have this interface slot. Refer to “[Interfaces](#)” on page 196 for details.

p/n	Description
G1369B or G1369-60002	Interface board (LAN)
G1369C or G1369-60012	Interface board (LAN)

NOTE

One board is required per Agilent 1260 Infinity instrument. It is recommended to add the LAN board to the detector with highest data rate.

NOTE

For the configuration of the G1369 LAN Communication Interface card refer to its documentation.

The following cards can be used with the Agilent 1260 Infinity modules.

Table 28 LAN Boards

Type	Vendor	Supported networks
Interface board (LAN) (p/n G1369B or G1369-60002) or Interface board (LAN) (p/n G1369C or G1369-60012)	Agilent Technologies	Fast Ethernet, Ethernet/802.3, RJ-45 (10/100Base-TX) <i>recommended for re-ordering</i>
LAN Communication Interface board (p/n G1369A or G1369-60001)	Agilent Technologies	Fast Ethernet, Ethernet/802.3, RJ-45 (10/100Base-TX) <i>(obsolete)</i>
J4106A ¹	Hewlett Packard	Ethernet/802.3, RJ-45 (10Base-T)

Table 28 LAN Boards

Type	Vendor	Supported networks
J4105A ¹	Hewlett Packard	Token Ring/802.5, DB9, RJ-45 (10Base-T)
J4100A ¹	Hewlett Packard	Fast Ethernet, Ethernet/802.3, RJ-45 (10/100Base-TX) + BNC (10Base2)

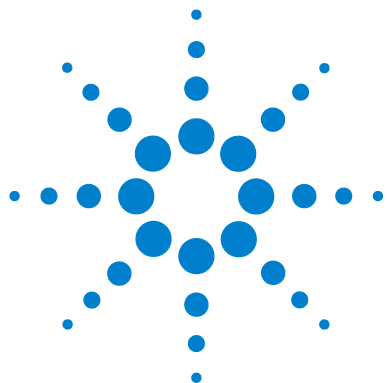
¹ These cards may be no longer orderable. Minimum firmware of these Hewlett Packard JetDirect cards is A.05.05.

Recommended LAN Cables

p/n	Description
5023-0203	Cross-over network cable, shielded, 3 m (for point to point connection)
5023-0202	Twisted pair network cable, shielded, 7 m (for point to point connection)

12 Hardware Information

Optional Interface Boards



13 Appendix

General Safety Information	212
General Safety Information	212
Safety Standards	212
Operation	212
Safety Symbols	214
The Waste Electrical and Electronic Equipment Directive	215
Lithium Batteries Information	216
Radio Interference	217
Sound Emission	218
Agilent Technologies on Internet	219

This chapter provides addition information on safety, legal and web.



General Safety Information

General Safety Information

The following general safety precautions must be observed during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions or with specific warnings elsewhere in this manual violates safety standards of design, manufacture, and intended use of the instrument. Agilent Technologies assumes no liability for the customer's failure to comply with these requirements.

WARNING

Ensure the proper usage of the equipment.

The protection provided by the equipment may be impaired.

→ The operator of this instrument is advised to use the equipment in a manner as specified in this manual.

Safety Standards

This is a Safety Class I instrument (provided with terminal for protective earthing) and has been manufactured and tested according to international safety standards.

Operation

Before applying power, comply with the installation section. Additionally the following must be observed.

Do not remove instrument covers when operating. Before the instrument is switched on, all protective earth terminals, extension cords, auto-transformers, and devices connected to it must be connected to a protective earth via a ground socket. Any interruption of the protective earth grounding will cause a potential shock hazard that could result in serious personal injury. Whenever it is likely that the protection has been impaired,

the instrument must be made inoperative and be secured against any intended operation.

Make sure that only fuses with the required rated current and of the specified type (normal blow, time delay, and so on) are used for replacement. The use of repaired fuses and the short-circuiting of fuse holders must be avoided.

Some adjustments described in the manual, are made with power supplied to the instrument, and protective covers removed. Energy available at many points may, if contacted, result in personal injury.

Any adjustment, maintenance, and repair of the opened instrument under voltage should be avoided whenever possible. When inevitable, this has to be carried out by a skilled person who is aware of the hazard involved. Do not attempt internal service or adjustment unless another person, capable of rendering first aid and resuscitation, is present. Do not replace components with power cable connected.

Do not operate the instrument in the presence of flammable gases or fumes. Operation of any electrical instrument in such an environment constitutes a definite safety hazard.

Do not install substitute parts or make any unauthorized modification to the instrument.

Capacitors inside the instrument may still be charged, even though the instrument has been disconnected from its source of supply. Dangerous voltages, capable of causing serious personal injury, are present in this instrument. Use extreme caution when handling, testing and adjusting.

When working with solvents, observe appropriate safety procedures (for example, goggles, safety gloves and protective clothing) as described in the material handling and safety data sheet by the solvent vendor, especially when toxic or hazardous solvents are used.

Safety Symbols

Table 29 Safety Symbols

Symbol	Description
	The apparatus is marked with this symbol when the user should refer to the instruction manual in order to protect risk of harm to the operator and to protect the apparatus against damage.
	Indicates dangerous voltages.
	Indicates a protected ground terminal.
	Indicates eye damage may result from directly viewing the light produced by the deuterium lamp used in this product.
	The apparatus is marked with this symbol when hot surfaces are available and the user should not touch it when heated up.

WARNING

A WARNING

alerts you to situations that could cause physical injury or death.

- Do not proceed beyond a warning until you have fully understood and met the indicated conditions.

CAUTION

A CAUTION

alerts you to situations that could cause loss of data, or damage of equipment.

- Do not proceed beyond a caution until you have fully understood and met the indicated conditions.

The Waste Electrical and Electronic Equipment Directive

Abstract

The Waste Electrical and Electronic Equipment (WEEE) Directive (2002/96/EC), adopted by EU Commission on 13 February 2003, is introducing producer responsibility on all electric and electronic appliances starting with 13 August 2005.

NOTE

This product complies with the WEEE Directive (2002/96/EC) marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.

Product Category:

With reference to the equipment types in the WEEE Directive Annex I, this product is classed as a Monitoring and Control Instrumentation product.



NOTE

Do not dispose off in domestic household waste

To return unwanted products, contact your local Agilent office, or see www.agilent.com for more information.

Lithium Batteries Information

WARNING

Lithium batteries may not be disposed-off into the domestic waste. Transportation of discharged Lithium batteries through carriers regulated by IATA/ICAO, ADR, RID, IMDG is not allowed.

Danger of explosion if battery is incorrectly replaced.

- Discharged Lithium batteries shall be disposed off locally according to national waste disposal regulations for batteries.
- Replace only with the same or equivalent type recommended by the equipment manufacturer.



WARNING

Lithiumbatteri - Eksplosionsfare ved fejlagtig håndtering.

Udskiftning må kun ske med batteri af samme fabrikat og type.

- Lever det brugte batteri tilbage til leverandøren.

WARNING

Lithiumbatteri - Eksplosionsfare.

Ved udskiftning benyttes kun batteri som anbefalt av apparatfabrikanten.

- Brukt batteri returneres apparatleverandøren.

NOTE

Bij dit apparaat zijn batterijen geleverd. Wanneer deze leeg zijn, moet u ze niet weggooien maar inleveren als KCA.

Radio Interference

Cables supplied by Agilent Technologies are screened to provide optimized protection against radio interference. All cables are in compliance with safety or EMC regulations.

Test and Measurement

If test and measurement equipment is operated with unscreened cables, or used for measurements on open set-ups, the user has to assure that under operating conditions the radio interference limits are still met within the premises.

Sound Emission

Manufacturer's Declaration

This statement is provided to comply with the requirements of the German Sound Emission Directive of 18 January 1991.

This product has a sound pressure emission (at the operator position) < 70 dB.

- Sound Pressure $L_p < 70$ dB (A)
- At Operator Position
- Normal Operation
- According to ISO 7779:1988/EN 27779/1991 (Type Test)

Agilent Technologies on Internet

For the latest information on products and services visit our worldwide web site on the Internet at:

<http://www.agilent.com>

Select Products/Chemical Analysis

It will provide also the latest firmware of the modules for download.

Index

8

8-bit configuration switch
without On-Board LAN 202

A

adapter required 111
Agilent Diagnostic software 95
Agilent Lab Advisor software 95
Agilent Lab Advisor 95
Agilent
on internet 219
ambient non-operating temperature 19
ambient operating temperature 19
analog signal 200
analog
cable 182
apg remote 200

B

battery
safety information 216
BCD board
external contacts 206
BCD
cable 187
bench space 18
bio-inert 128
materials 10
board
HP JetDirect card 208
boards
LAN card 208

C

cable
analog 182
BCD 187
CAN 189
external contact 190
LAN 189
remote 184
RS-232 191
cables
analog 180
BCD 180
CAN 181
external contact 181
LAN 181
overview 180
remote 180
RS-232 181
calib delay vol two peaks 110
CAN
cable 189
cleaning 129
cluster partner lost during analysis 114
Communication settings
RS-232C 203
compensation sensor open 103
compensation sensor short 104
condensation 18
could not find a valid next position 115
D
defect on arrival 24
Diagnostic software 95
dimensions 19

E

electrical connections
descriptions of 194
electronic waste 215
electrostatic discharge (ESD) 128, 160
EMF
early maintenance feedback 12
error messages
calib delay vol two peaks 110
cluster partner lost during
analysis 114
could not find a valid next
position 115
lost CAN partner 102
motor temperature 113
remote timeout 101
vessel stuck to needle 114
error messages
adapter required 111
arm movement 107
compensation sensor open 103
compensation sensor short 104
fan failed 104
ignition without cover 105, 105
leak sensor open 103
leak sensor short 102
leak 105
shutdown 101
timeout 100
external contact
cable 190
external contacts
BCD board 206

F

fan failed 104
 firmware
 updates 162, 162
 upgrade/downgrade 162
 upgrade/downgrade 162
 frequency range 19

G

general error messages 100

H

HP JetDirect card 208
 humidity 19

I

installation
 bench space 18
 instrument layout 13
 interfaces 196
 internet 219

L

LAN
 cable 189
 communication interface board 208
 leak sensor open 103
 leak sensor short 102
 leak 105
 line frequency 19
 line voltage 19
 lithium batteries 216
 lost CAN partner 102

M

maintenance
 definition of 126

feedback 12
 replacing firmware 162, 162

materials

bio-inert 10

message

ignition without cover 105, 105
 remote timeout 101

N

non-operating altitude 19
 non-operating temperature 19

O

operating Altitude 19
 operating temperature 19

P

packaging
 damaged 24
 physical specifications 19
 power consumption 19
 power supply indicator 93
 power consideration 16
 power cords 17

R

radio interference 217
 remote
 cable 184
 repairs
 replacing firmware 162, 162
 RS-232C
 cable 191
 communication settings 203

S

safety class I 212
 safety information

lithium batteries 216

safety

general information 212
 standards 19
 symbols 214

shutdown 101

site requirements

power cords 17
 solvent information 85

sound emission 218

special interfaces 201

special settings

boot-resident 205
 forced cold start 205

specification

physical 19

status indicator 94

T

temperature sensor 105
 timeout 100
 troubleshooting
 error messages 99, 92
 status indicators 92, 93

V

voltage range 19

W**waste**

electrical and electronic
 equipment 215

WEEE directive 215

weight 19

In This Book

This manual contains technical reference information about the Agilent 1260 Infinity Bio-inert Analytical-scale Fraction Collector G5664A.

The manual describes the following:

- introduction to the fraction collector,
- site requirements and specifications,
- installing the fraction collector,
- using and optimizing,
- troubleshooting and test functions,
- maintenance,
- parts and materials,
- hardware,
- cables, and
- safety information.

© Agilent Technologies 2011-2012

Printed in Germany
05/2012



G5664-90001 Rev. B



Agilent Technologies