

Agilent 1220 Infinity II LC Mobile Upgrade Kit

Installation Note

This note describes the procedures to install an Mobile Upgrade kit to an existing Agilent 1220 Infinity II LC.

Contents

Introduction	2
Application Examples	2
Site Requirements and Specifications	3
Site Requirements	3
Specifications	6
Installation	7
Cautions and Warnings	7
Unpacking the Module	7
Installing the Handle Plate to the 1220 Infinity II LC Instrument	10
Installing the Solvent Selection Valve Support	12
Installing the Mobile Solvent Compartment	13
Installing the Column Wire Mesh	15
Installing the Attenuation Unit to the Bench	15
Securing the LC to the attenuation unit	17
General Operation Instructions	19
Consumables	20



Introduction

The Agilent 1220 Infinity II Mobile LC Solution is an Agilent 1220 Infinity II LC with integrated variable wavelength or diode array detector and a mobile upgrade kit that enables transportation. The 1220 Infinity II LC Mobile Upgrade Kit consists of functional parts that allows the Agilent 1220 Infinity II LC to be mounted into a mobile lab so it can be moved to different locations to access remote measurement sites. The main component is the attenuation unit which acts as a shock absorber to protect the instrument during transit or from influences of operators moving in the mobile lab. A solvent bottle unit keeps the bottles fixed to the instrument. A column wire mesh keeps the column safe in the column heater and the 'mobile solvent compartment' secures the solvent bottles during operation of the instrument.

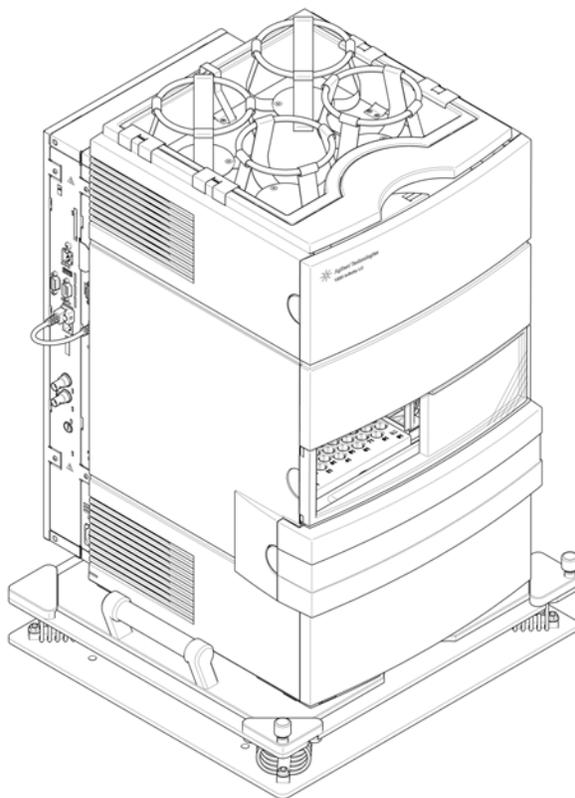


Figure 1 Agilent 1220 Infinity II Instrument with attached mobile kit

Application Examples

When samples of interest are locally remote from the lab, the lab has to come to the samples. GC and GC/MS based mobile labs have been around for quite some time but the need for mobile LCs has become a more burning issue recently. Mobile LC applications range from analysis of non-volatile pesticides in environmental and food applications, monitoring of drugs or testing of counterfeit drugs. It can also be used in teaching, when a mobile lab goes to students residing in remote areas. The diode array detector integrated in the Agilent 1220 Infinity II LC enables spectra analysis for purity identification and comparison with self-created library data.

Site Requirements and Specifications

Site Requirements

Bench Space

The attenuation unit has a footprint of 480 mm (18.9 in) in width and 445 mm (17.5 in) in depth.

The overall room specified for operation (width x depth x height) of 560 mm (22.1 in) x 665 mm (26.2 in) x 970 mm (38.2 in) already includes the space for a complete 1220 Infinity LC Instrument mounted onto the attenuation unit and considering the necessary room for air circulation, electrical connections and safety area for the possible movement of the instrument on the attenuation unit.

For a comfortable handling of the solvent bottles additionally 20 cm (7.9 in) should be kept free above the bottles.

See [Figure 2](#) on page 3 for more details.

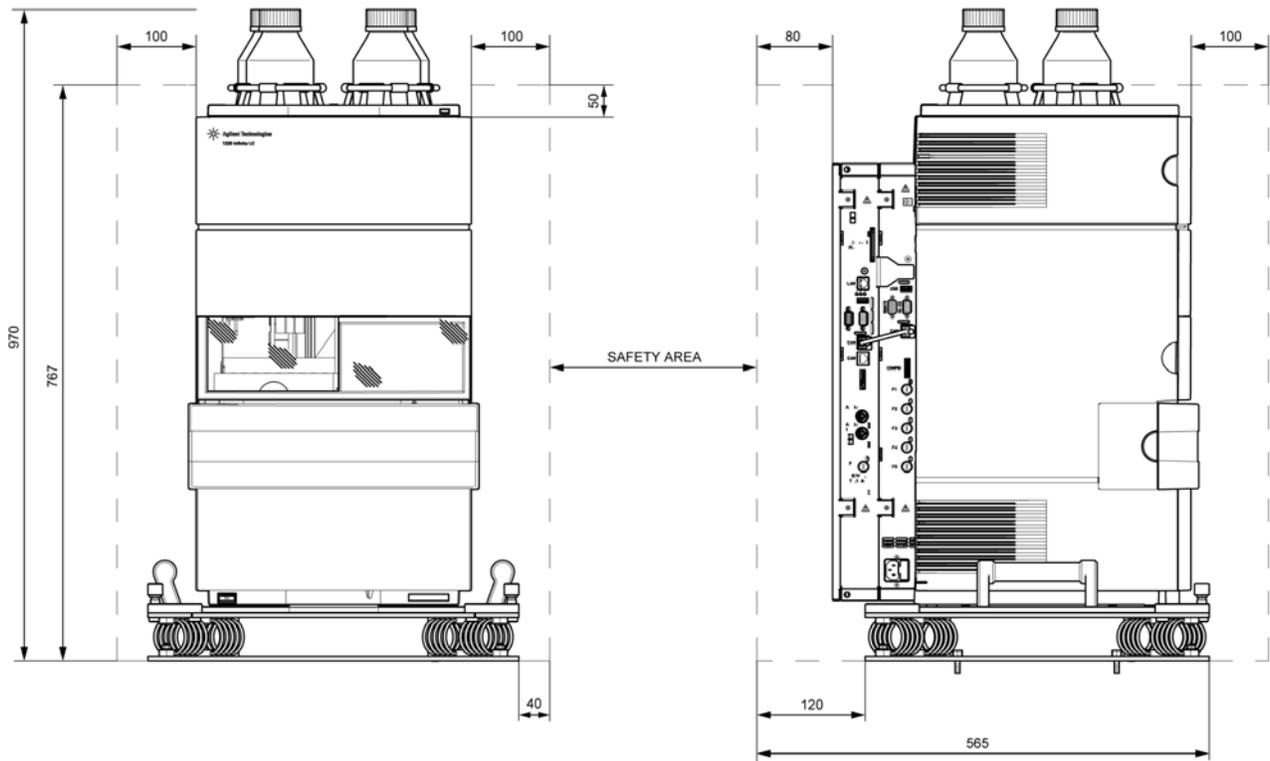


Figure 2 Required space for a 1220 Infinity II LC Instrument installed to a mobile upgrade kit (numbers given in mm)

Bench Material and Strength

A proper bench must be chosen under the aspect that it must bear the combined weight of the instrument with solvent bottles and the mobile kit.

The bench material must be suitable for a proper fastening solution.

Fastening Solution

Because of the expected varieties in bench material and strength the fastening material is not included in the kit.

A suitable fastening solution (screws or bolts) must be chosen to secure the attenuation unit to the bench of the mobile lab.

NOTE

This solution must be strong enough to withstand the forces that will result from the possible accelerations that can occur in a moving vehicle (app. 2000 N per screw, based on a calculation with 1 G acceleration and the weight of a G4294B).

NOTE

The head of the chosen screws or bolts must not stand out more than 15 mm on top of the attenuation unit's bottom plate.

The attenuation unit comes with four 8.5mm (0.33 in) diameter holes prepared in the lower plate and four 20 mm (0.79 in) diameter holes located directly above them in the top plate. This allows the use of a 14 mm (9/16 in) nut for tightening the screws or bolts.

See [Figure 3](#) on page 5 for the layout information of the holes that must be prepared to the bench.

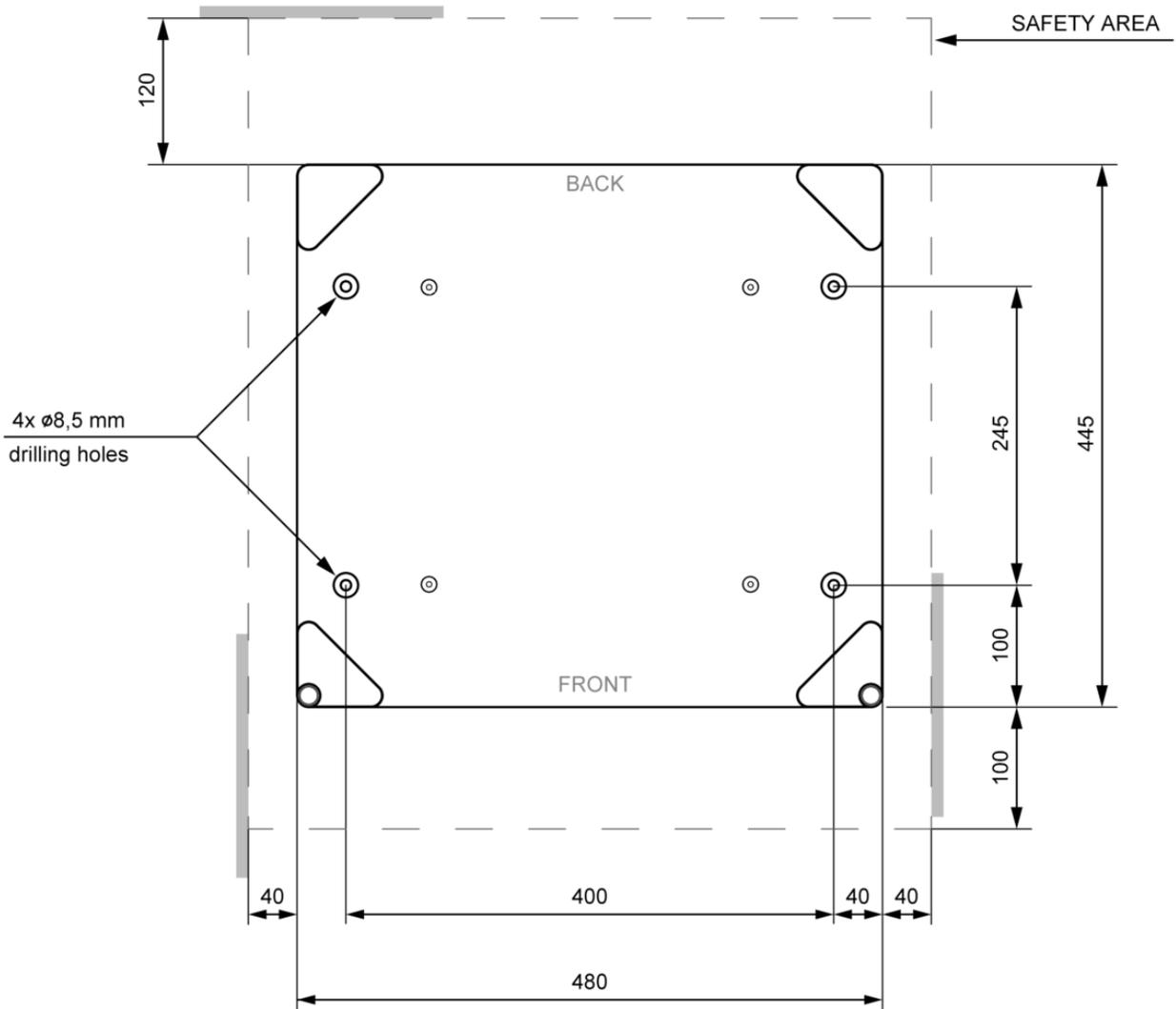


Figure 3 Dimensions of the attenuation unit

Specifications

Physical Specifications

Table 1 Physical Specifications

Type	Specification	Comments
Weight (all parts)	14.2 kg (31.3 lbs)	
Dimensions of the attenuation unit (height x width x depth)	73 x 480 x 445 mm (2.9 x 18.9 x 17.5 in)	See “Installing the Attenuation Unit to the Bench” on page 15 for more information regarding safety space.

Performance Specifications

The performance specifications show details about the performed test patterns.

Table 2 Performance Specifications

Type	Specification	Comments
Vibration Test	Outdoor	Agilent internal test guidelines

Installation

Cautions and Warnings

WARNING

Personal injury or damage to the system

The system is bulky and heavy.

→ Lift the instrument with at least two persons.

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.

CAUTION

Damage to the system

If the handle plate is not fixed precisely, the mobile system may tumble or slide off the bench.

→ Use the handle plate only if it is fixed correctly to the instrument.

NOTE

Please note that the design of the *1220 Infinity II mobile solution* is intended for protecting the instrument from influences during transit. Analytical measurement however has to be conducted under stationary conditions.

NOTE

Operating the 1220 Infinity II LC in a mobile lab environment can require special equipment or facilities to comply with local safety regulations (e.g. special solvent waste and storage containers). Please verify that these regulations are met before you operate the 1220 Infinity II LC in your mobile lab.

Unpacking the Module

Damage to the module

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 mobile upgrade kit. The delivery checklist is shown below. Please report missing or damaged parts to your local Agilent Technologies sales and service office.

p/n	Description
G4292-60001	Mobile solvent compartment Includes 4 prefixed bottle holders and 4 Leak Tub brackets to fix it to the module's solvent compartment
G4292-60000	Attenuation unit Shock absorbing unit to mount on the Mobile Lab's bench
G4292-60003	Handling plate Pre-connected to the Attenuation unit (see " Installing the Handle Plate to the 1220 Infinity II LC Instrument " on page 10 for further instructions)
G4292-40000	Column wire mesh To protect the column
G4292-00001	SSV support Needs to be installed when a solvent selection valve is installed

NOTE

The required fastening bolts or screws are not delivered with the mobile upgrade kit and a proper solution must be chosen. See for more details.

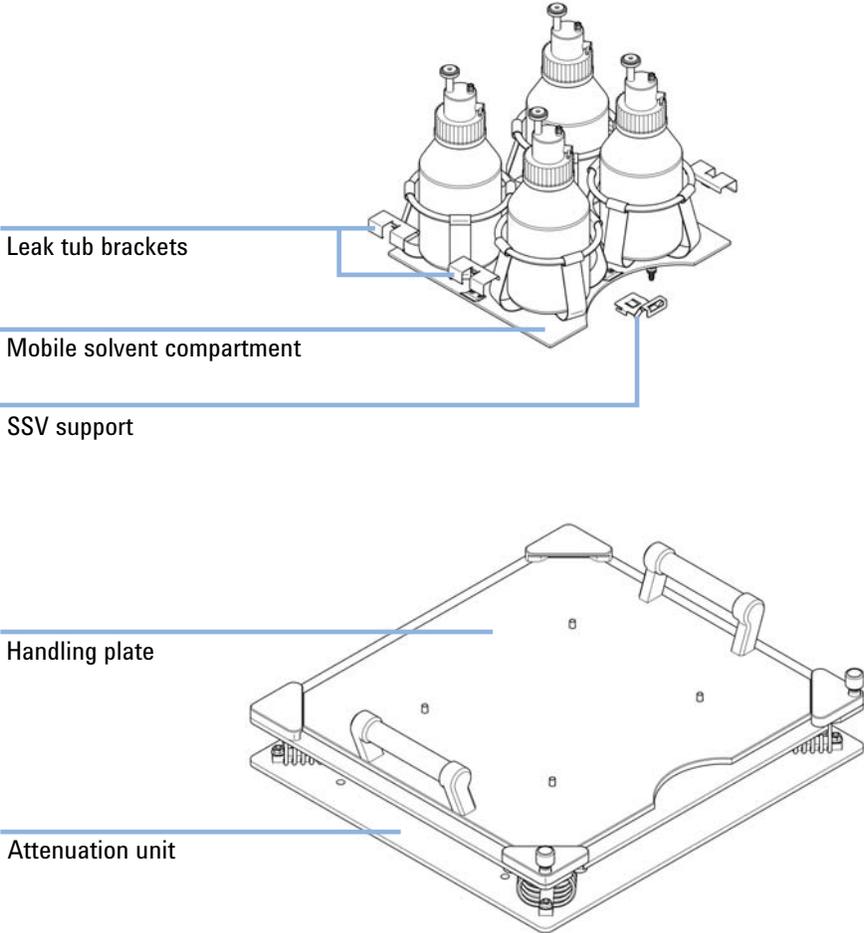


Figure 4 Mobile Kit Overview

Installing the Handle Plate to the 1220 Infinity II LC Instrument

WARNING

Personal injury or damage to the system

The system is bulky and heavy.

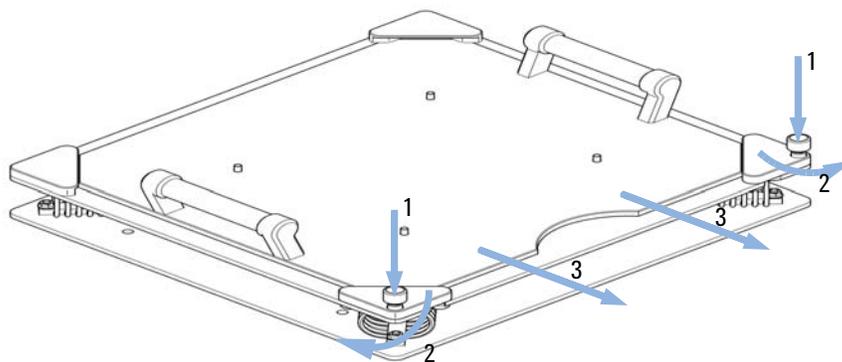
→ Lift the instrument with at least two persons.

1 Lay the 1220 Infinity II LC Module carefully on its rear plate.

NOTE

For the G4290B it is necessary either to position the bottom of the instrument slightly free over the bench's edge or to lay it onto the carton box of the solvent compartment.

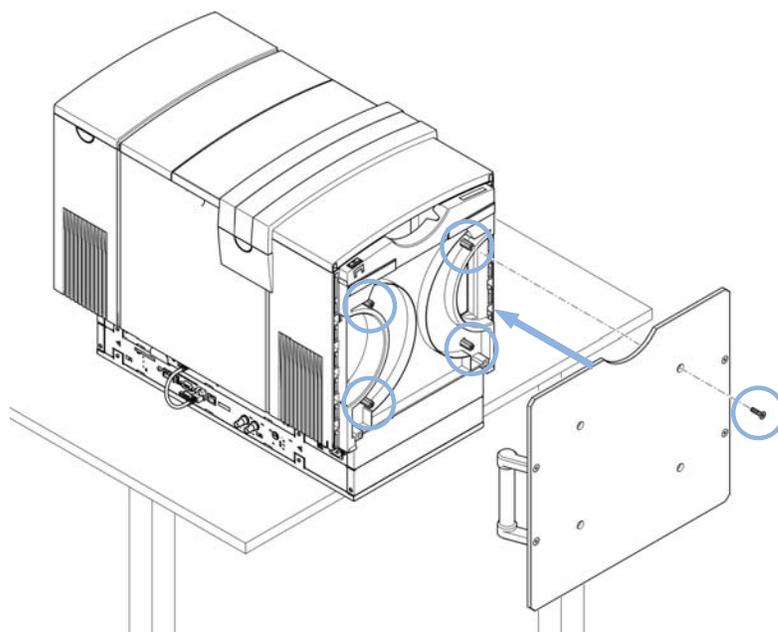
2 Push on the two lock pins (1) to release the front fixing angles (2) and separate the handling plate (3) from the rest of the attenuation unit.



3 Use the four delivered screws to mount the handle plate to the bottom of the 1220 Infinity II LC instrument. The holes for the screws are already prepared in the bottom of the instrument.

NOTE

Pay attention that the side with the concave cut-out section is orientated to the instrument front.



- 4 Verify that the screws are fixed and carefully bring the instrument back into the upright position.

Installing the Solvent Selection Valve Support

When The Solvent Selection Valve Support needs to be installed if a solvent selection valve (SSV) is used with the instrument.

Parts required	p/n	Description
	G4292-00001	SSV support (delivered as part of the G4292-60000 attenuation unit)

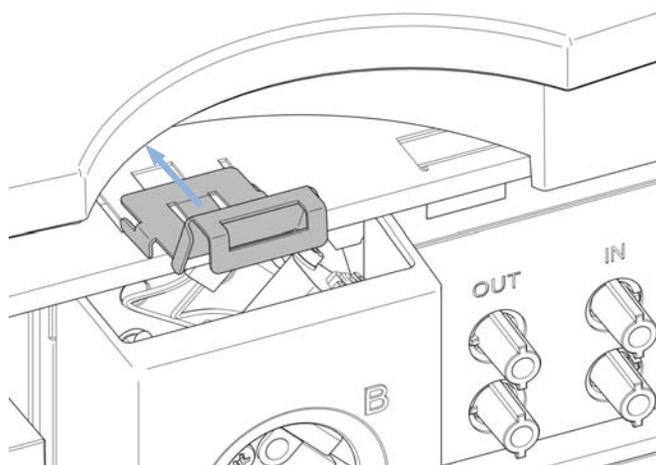
Preparations If SSV is already installed:

- Unplug the SSV.
- Clear the solvent compartment.
- Remove the whole top cover with the SSV still attached to it.

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.

NOTE When the SSV support is installed, the SSV cannot get attached to the top cover without removing the top cover.

- 1 Take the SSV-Support and slide it into the slits in the metal cover above the gradient valve.



- 2 If the SSV was already installed:
 - Reinstall the top cover.
 - Connect the SSV.OR
If the SSV was not already installed:
 - Remove the top cover.
 - Fix the SSV to the top cover.
 - Reinstall the top cover.
 - Connect the SSV.

Installing the Mobile Solvent Compartment

Parts required	p/n	Description
	G4292-60001	Mobile solvent compartment including 4 special brackets

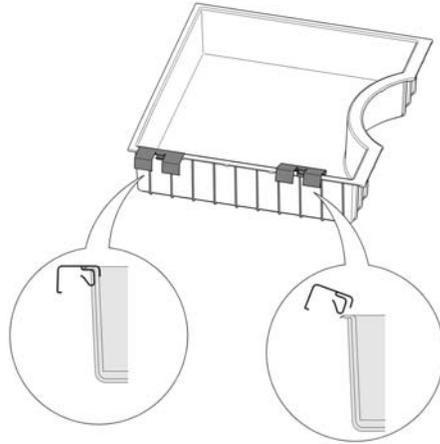
WARNING

Sharp metal edges

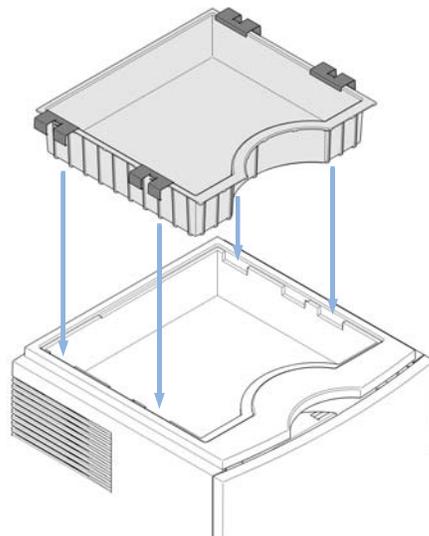
Sharp-edged parts of the equipment may cause injuries.

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

- 1 Take out the grey solvent bottle container of the instrument's top cover and attach the 4 leak tub brackets to it (two on each side). Look at the instrument's top and identify the referring spaces (where the arrows in the picture of step 2 on page 13 point to). This is the dedicated position for middle part of the clamps. This information is necessary for a correct positioning of the clamps. It is hard to correct their position once they are installed.



- 2 Now take the leak tub and reinstall it to the top cover. Verify that the leak tub brackets properly hook in underneath the top cover on the instrument's side.

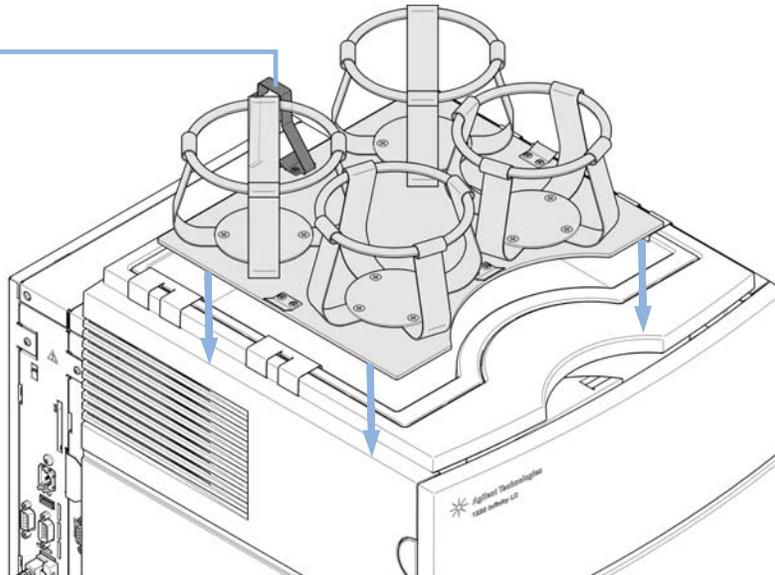


Installation

Installing the Mobile Solvent Compartment

- 3 Take the mobile solvent compartment and push it into the solvent container. The two metal clips will grip into the solvent container's plastic. On the rear of the mobile solvent compartment is the grounding clamp which needs to hook into the metal sheet between the board extension and the main assembly.

Grounding clamp



Installing the Column Wire Mesh

Parts required	p/n	Description
	G4292-40000	Column wire mesh

- 1 Put the column wire mesh with its open side facing to the top into the column heater assembly.
- 2 Position the column in the Column wire mesh so that it is fixed properly.

Installing the Attenuation Unit to the Bench

The bench holding the instrument requires 4 screws or bolts to secure the attenuation unit of the 1220 Infinity II Mobile LC Kit to the bench. This fastening material is not provided from Agilent because of the unpredictable situation of bench material and dimension.

Tools required	Description
	Wrench (size depending on chosen fastening material)

Parts required	p/n	Description
	G4292-60000	Attenuation unit

Preparations

Prepare the appropriate fastening material (4 proper screws or bolts) and corresponding tools for attaching the attenuation unit onto the desired surface.

Four 8.5 mm (0.33 in) holes are prepared on the lower plate of the attenuation unit and four equivalent 20 mm (0.79 in) holes in the top plate above allowing the use of a 14mm nut.

The screws or bolts must withstand a maximum drag force of 2000 N and their head should not stand out more than 15 mm out of the lower plate.

See figure in step 2 on page 16 for the layout location of these holes that must be provided in the bench top.

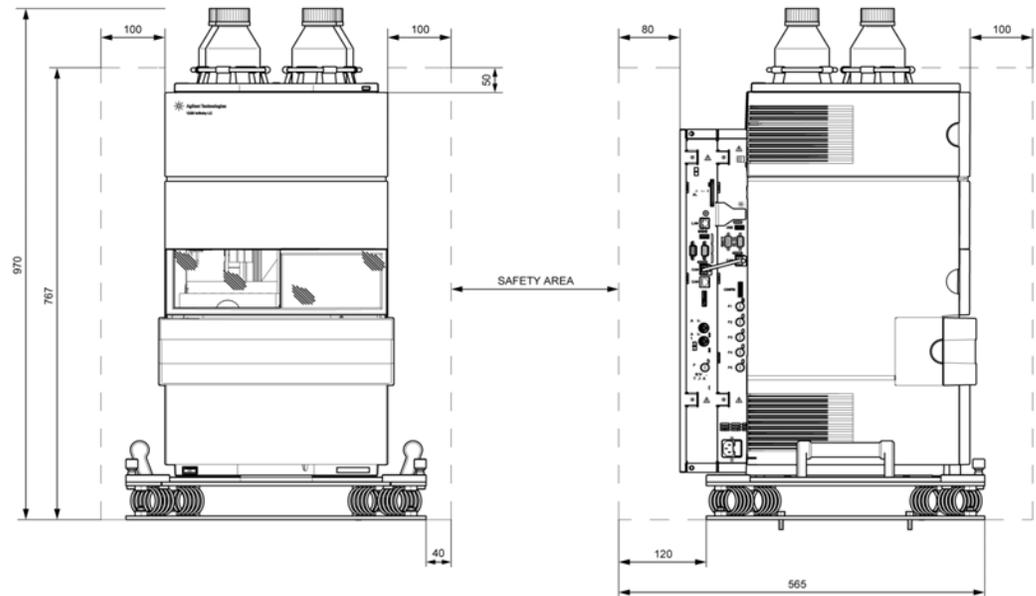
NOTE

The fastening material is not part of this kit.

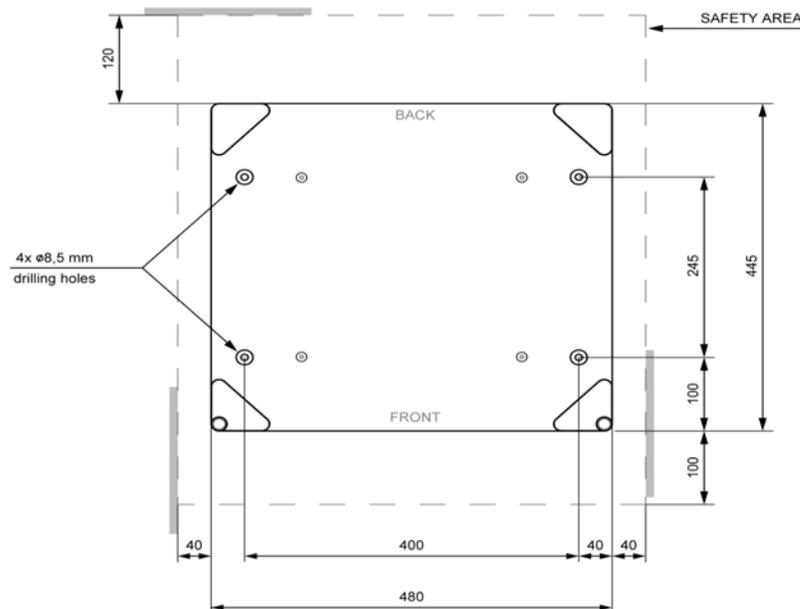
Installation

Installing the Attenuation Unit to the Bench

- 1 Identify the final position of the instrument on the bench according to the specifications given below.



- 2 Drill the 4 holes in the bench. Either use the attenuation unit as a model or prepare the holes according to the figure below.



- 3 Position the attenuation unit on the bench. The snap fasteners must be oriented to the front and the prepared holes in the attenuation unit must stand above the drill-holes in the bench.
- 4 Use the chosen screws or bolts to fix the attenuation unit to the bench.

Securing the LC to the attenuation unit

- When** Whenever the instrument is installed onto the attenuation unit in the mobile lab.
- Preparations** Install the handle plate to the 1220 Infinity II LC Instrument (see [“Installing the Handle Plate to the 1220 Infinity II LC Instrument”](#) on page 10).
Install the attenuation unit to the bench (see [“Installing the Attenuation Unit to the Bench”](#) on page 15).

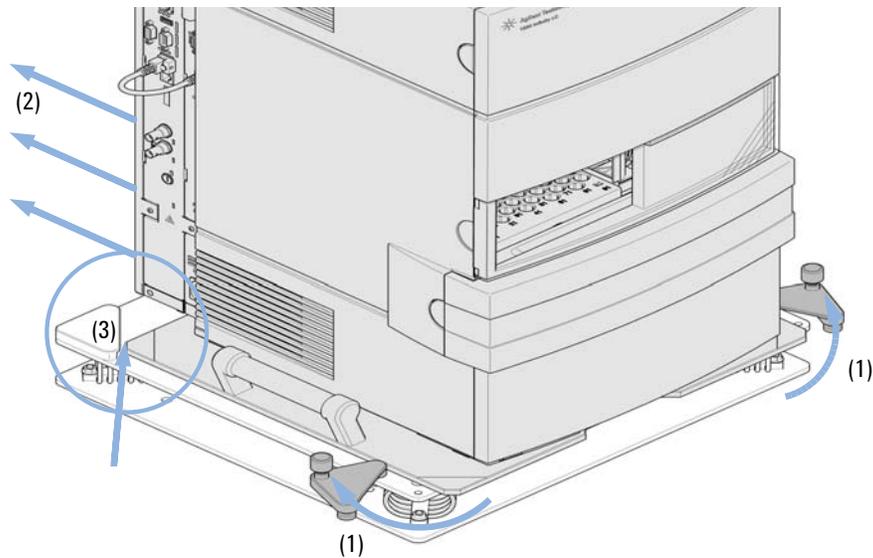
WARNING

Personal injury or damage to the system

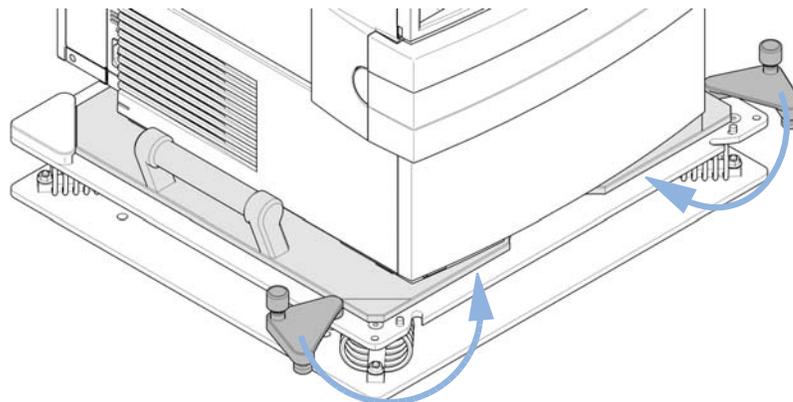
The system is bulky and heavy.

→ Lift the instrument with at least two persons.

- 1 Open the fixing angles (1) and slide the instrument on the handling plate onto the attenuation unit (2). Verify that it is completely fixed by the two rear retaining brackets (3).



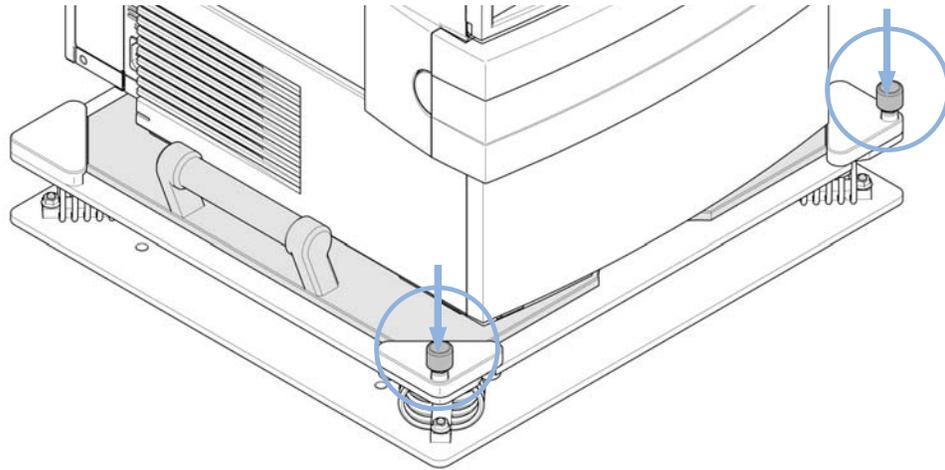
- 2 Close the two front retaining brackets.



Installation

Securing the LC to the attenuation unit

- 3 Secure the front retaining brackets by pressing the arrester button.



NOTE

The two front retaining brackets must always get closed and secured by pressing the arrester button. They will not secure the instrument if not properly closed.

General Operation Instructions

NOTE

Please note that the design of the *1220 Infinity II mobile solution* is intended for protecting the instrument from influences during transit. Analytical measurement however has to be conducted under stationary conditions.

NOTE

Operating the 1220 Infinity II LC in a mobile lab environment can require special equipment or facilities to comply with local safety regulations (e.g. special solvent waste and storage containers). Please verify that these regulations are met before you operate the 1220 Infinity II LC in your mobile lab.

Preparing transit and relocation

transit: 1220 Infinity II Instrument installed on the attenuation unit while the mobile lab is moving.

relocation: Moving the 1220 Infinity II Instrument to or from the mobile lab.

- Park the arm of the autosampler before relocation or transit.
- Remove the solvent bottles from the solvent cabinet and store them in a proper safety storage.
- Verify that the retaining brackets on the attenuation unit are properly locked when the instrument is on the attenuation unit (see “[Securing the LC to the attenuation unit](#)” on page 17).
- Store the sample vials in a safe place during relocation and transit. The sample tray is not a proper storage place.

General Operation

Use sealed solvent bottle caps and waste containers to prevent dangerous vaporization. More information can be found in the chapter “[Consumables](#)” on page 20 in this manual and also on the internet: www.agilent.com/chem/safetycaps

Consumables



Safety Cap I with 1 stopcock - GL45 (5043-0225)
(Comes with Venting valve (5043-0232))



Venting valve (5043-0232) with 1 µm PTFE membrane



10 L waste can S60 with Safety Cap IV 4 ports + 3 port collector (5043-0243)
Comes with a Safety Cap IV 4 ports - 1 leak port - S60 (5043-0227) including 4
ports and 1 leak port.
The Charcoal filter (48 g) (5043-0230) must be ordered seperately.



Charcoal filter (48 g) (5043-0230)



G4292-90001

Part Number: G4292-90001

Edition: 01/2017
Printed in Germany

© Agilent Technologies, Inc 2012-2016,
2017

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