

Agilent InfinityLab LC

Companion

Technical Note

This technical note describes the installation of the G7109A Agilent InfinityLab Companion Docking Station and Touchscreen Tablet (optional) for use with the G7108AA InfinityLab LC Companion.

Contents

Before You Start 2

Install the InfinityLab Companion Docking Station 3

Install the InfinityLab Companion Touchscreen Tablet 5

Install a non-Agilent Touchscreen Tablet 7

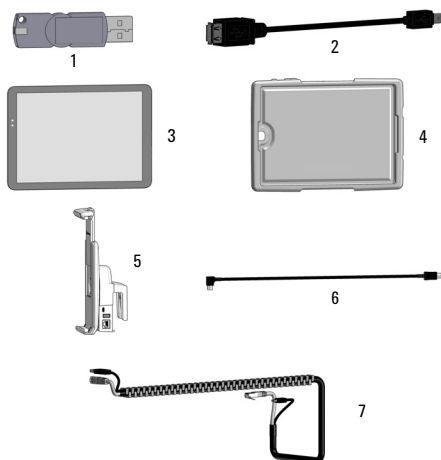
Electrical Interconnection 9



Agilent Technologies

Before You Start

Hardware required



- 1 InfinityLab Companion Docking Station (G7109A, Item 5), including Cable Set (Items 2, 6, and 7)
- 2 Infinity II Solvent Cabinet (sold separately, not shown)
- 3 Optional: InfinityLab Companion Touchscreen Tablet (G7109A#250, Item 3)
- 4 Optional: InfinityLab Companion Safety Case for (G7109A#252, Item 4)

Additionally, to enable the LC instrument to work with the InfinityLab Companion Touchscreen Tablet, the following item is required:

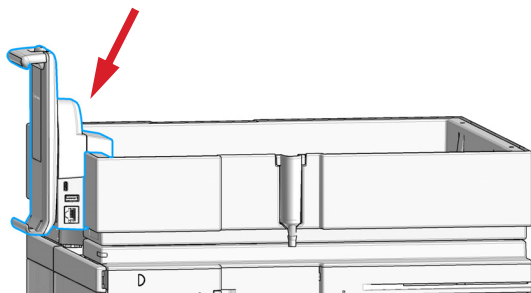
- InfinityLab LC Companion USB-Dongle (G7108AA, Item 1), including USB Adapter Cable

Preparations

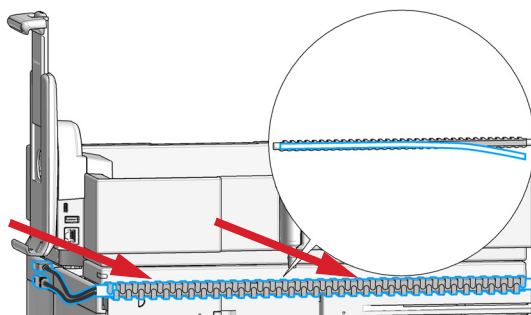
- Make sure that the InfinityLab LC Companion USB-Dongle (G7108AA, sold separately) is installed and activated for use on at least one of the modules of your LC instrument (see *Agilent InfinityLab Companion Software Installation Technical Note D0000195*).
- Make sure that the instrument on top of which the InfinityLab Docking Station will be installed, is turned off, before starting the installation.

Install the InfinityLab Companion Docking Station

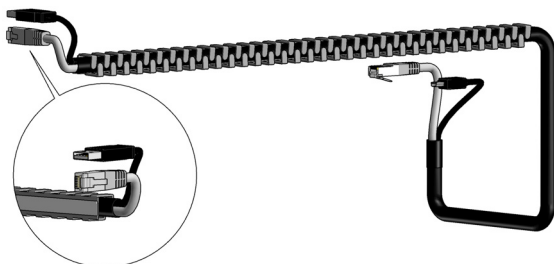
- 1 Attach the docking station to the solvent cabinet.



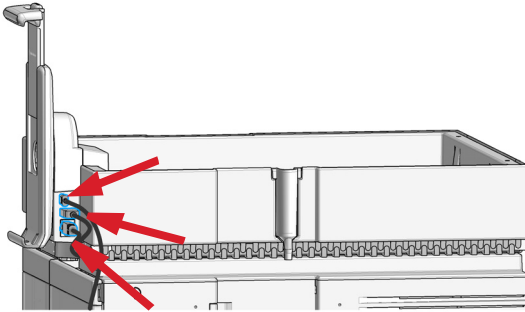
- 2 Fix the CAN/USB cable assembly centered with the adhesive tape along the right side of the solvent cabinet.



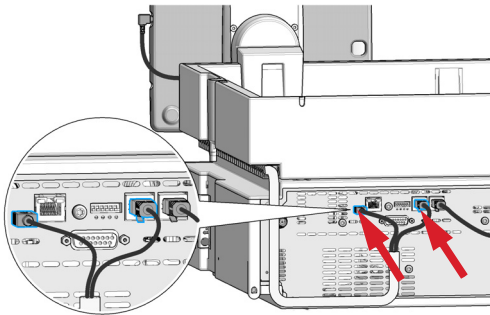
The USB A / CAN connector should be on the left side.



- 3** Plug the Micro-USB B power cable and the USB A/CAN cables into the ports at the docking station. If the cables are too long at the docking station, pull the cable assembly towards the rear through the channel.



- 4** Plug the CAN and the USB mini-b cables into the respective ports of the module holding the LAN connection and an activated InfinityLab LC Companion license.

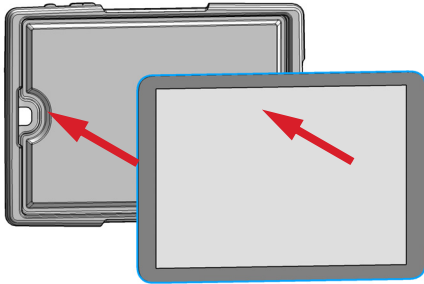


Install the InfinityLab Companion Touchscreen Tablet

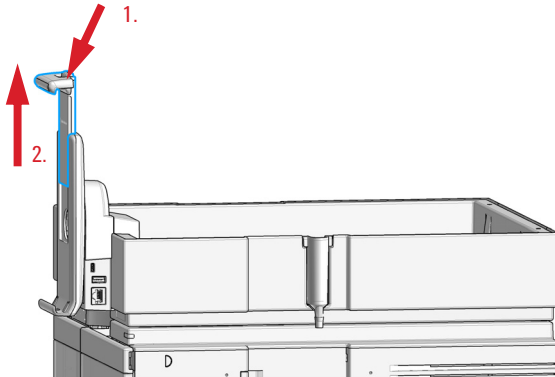
- 1 Insert the tablet into the Safety Case (Option #252, recommended)

OR

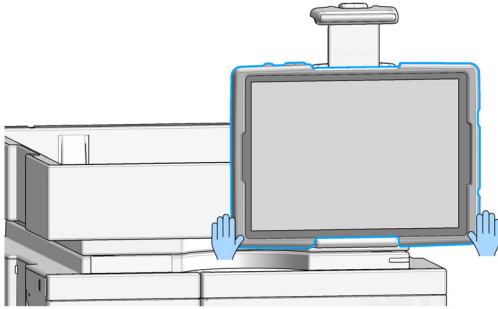
attach the rubber pads (included in G7109A) to the top and bottom of the tablet.



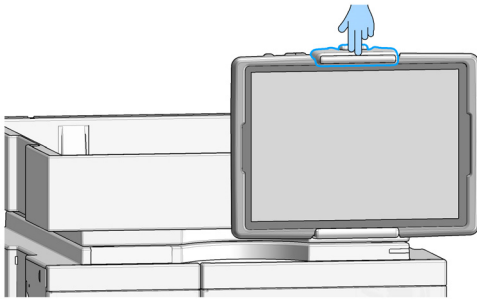
- 2 Press on the lid to release the slider (1.). Lift the slider up (2.).



3 Insert the tablet into the bottom of the docking station.

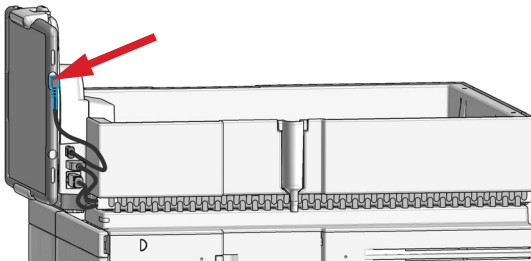


4 While pressing the release button, press the slider down to fix the tablet in the docking station.



5 Rotate the slider and tablet to achieve either portrait or landscape mode.

6 Connect the curved end of the micro-USB power connector to the tablet.



7 Power on the instrument.

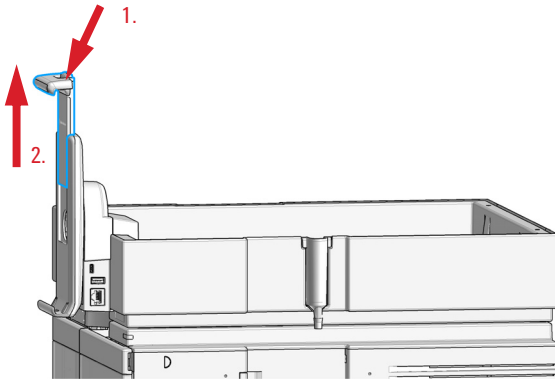
8 Power on the tablet.

Install a non-Agilent Touchscreen Tablet

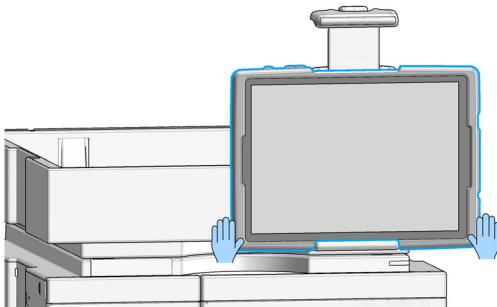
NOTE

When using a device other than the InfinityLab Companion Touchscreen Tablet, USB connectivity to the LC instrument is not supported. In this case, a connection to the LC instrument can only be achieved by using an existing WiFi connection.

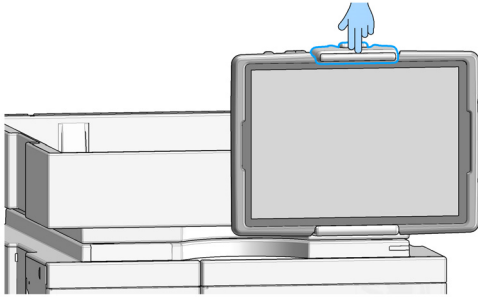
- 1 Attach the rubber pads included in G7109A InfinityLab Companion Docking Station to the top and bottom of your tablet device.
- 2 Press on the lid to release the slider (1.). Lift the slider up (2.).



- 3 Insert the tablet into the bottom of the docking station.



4 While pressing the release button, press the slider down to fix the tablet in the docking station.



5 Make sure that the tablet is attached firmly to the Docking Station.

Electrical Interconnection

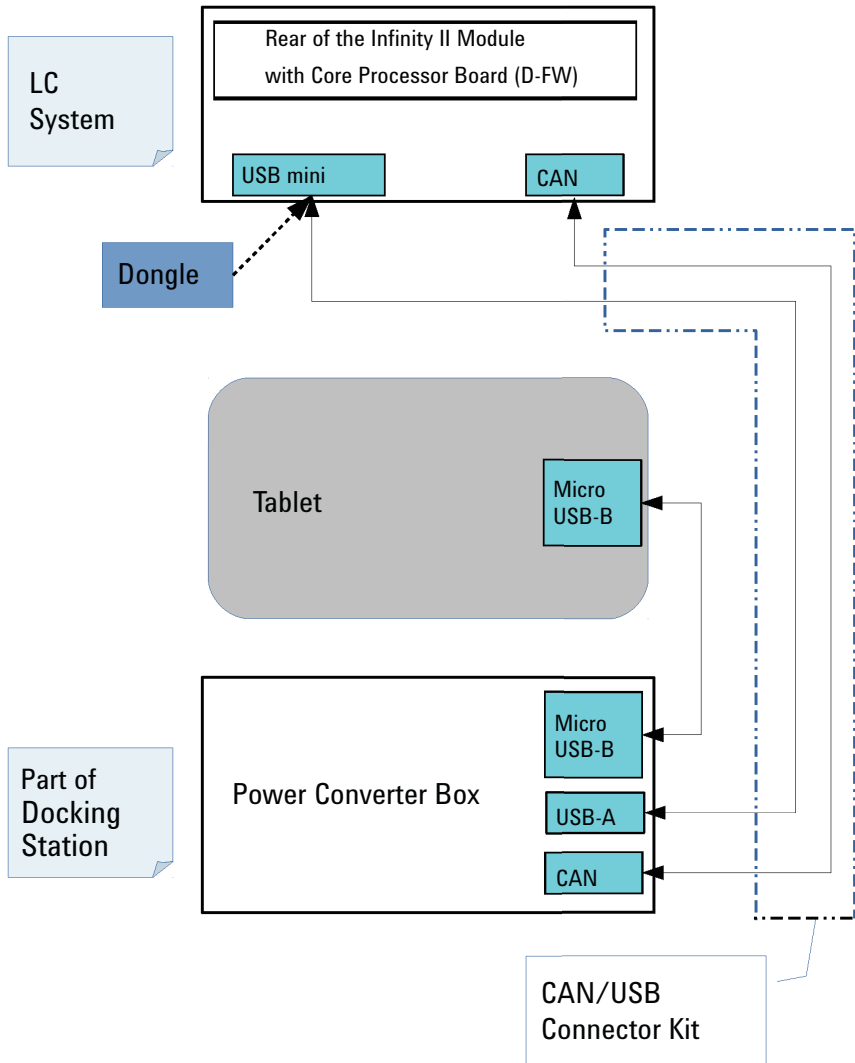


Figure 1 Electrical interconnection



G7108-90000

Part Number:
G7108-90000 Rev. D
Document number:
D0000216
Edition: 07/2019
Printed in Germany

© Agilent Technologies, Inc
2017-2019

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