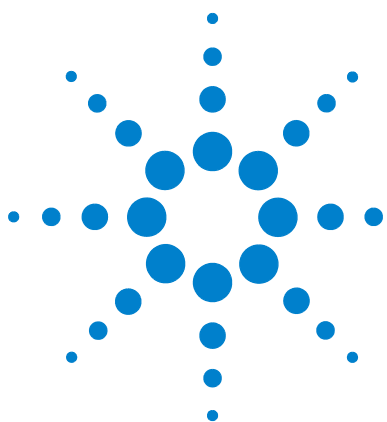




• Agilent 3100 OFFGEL Fractionator

Installation Quick Guide with Windows XP™

This instruction assist you to set up a connection from a PC with the operating system Windows XP professional to the Agilent 3100 OFFGEL Fractionator using a cross over LAN cable.



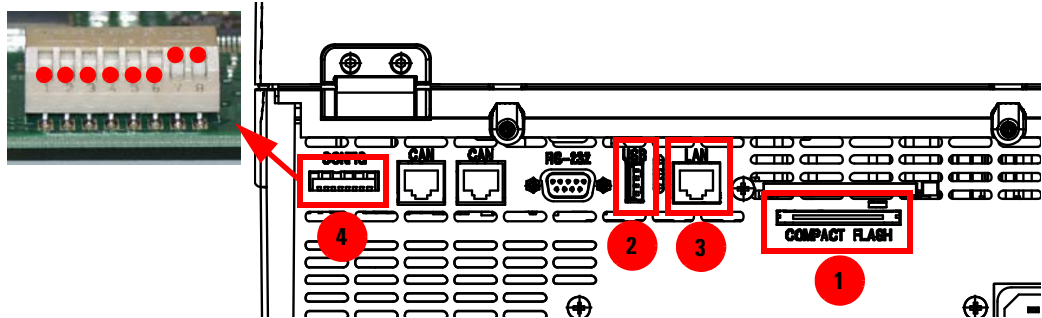
Tasks

- Preparation of the Agilent 3100 OFFGEL instrument.
- Assignment of an IP address to the PC.
- Access the instrument with Internet Explorer.



Agilent Technologies

Preparation of the Agilent 3100 OFFGEL Instrument



At the back of the instrument:

- ☐ Insert flash card to the instrument. (1)
- ☐ Insert USB stick. (2)
- ☐ Connect Cross-Over LAN cable. (3)
- ☐ Verify Configuration Switches: (4)
Switches N° 1 to 6 → down
Switches N° 7 and 8 → up
- ☐ Power the instrument on to apply this configuration.

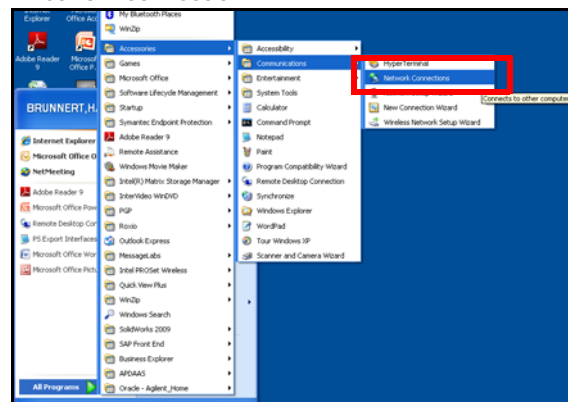
NOTE

This configuration enables the default IP address of the Agilent 3100 OFFGEL fractionator

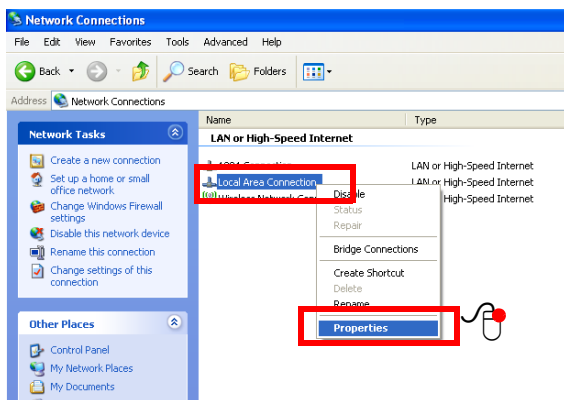
Assignment of a IP address to the PC

- 1 Plug Cross-Over LAN cable into the LAN connector of the PC.
- 2 Switch on the PC and Log on as administrator

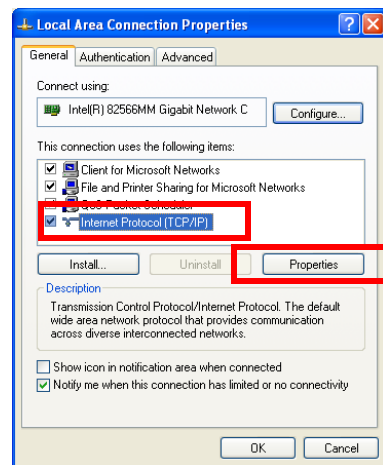
- 3 Select **All Programs > Accessories > Communications > Network connection**



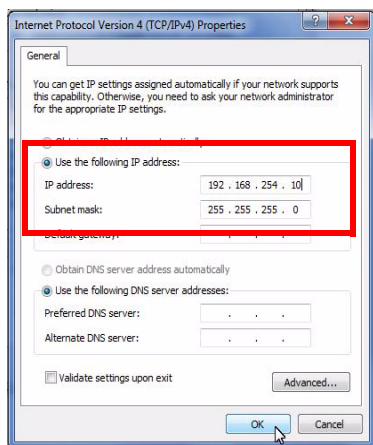
- 4 Select **Local Area Network**.
- 5 Click with the right mouse button
- 6 Select **Properties**



- 7 Select **Internet Protocol (TCP/IP)**.
- 8 Select **Use the following IP address**.
- 9 Click onto **Properties**.



- 10 Enter the following information:
IP address 192.168.254.10
Subnet mask 255.255.255.0
- 11 Select **Ok** to finish



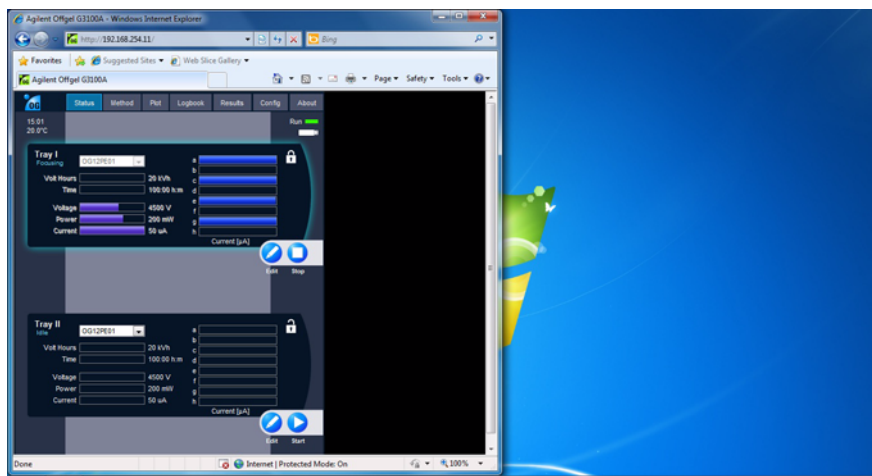
- 12 Close remaining windows.

NOTE

Your PC is now configured to operate within the same network address range as the Agilent 3100 OFFGEL fractionator.

Access the instrument with Internet Explorer

- 1 Start the Internet Explorer
- 2 Type IP address 192.168.254.11 into the address line.



NOTE

This address reflects the default IP address of the Agilent 3100 OFFGEL fractionator. It is different from the IP address of your PC.

Requirements and Specifications

The Instrument requires firmware B.06.30 or higher. Internet Explorer IE 7.0 and 8.0 support the web based user interface of the Agilent 3100 OFFGEL fractionator.

Network connection

To connect the Agilent 3100 OFFGEL fractionator to a local network request a static IP address from your administrator. Access the instrument first with the steps given above. Then set all supplied values under **Config** of the web-based User Interface, selecting **Set**. Power the instrument off and connect it to your local network with a standard LAN cable. Reconfigure the configuration switches to switch 7 up and all others down and restart the instrument. Your PC needs to be reconfigured to access the network and the instrument. For details refer to the 3100 OFFGEL user manual.

NOTE

Note down the initial Network configuration of the PC for your own reference.

