



Agilent G1369A LAN Interface



User Manual



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2003

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

G1369-90000

Edition

10/08/2003

Printed in Germany

Agilent Technologies, Deutschland GmbH
Hewlett-Packard-Strasse 8
76337 Waldbronn

Microsoft[®] is a U.S. registered trademark of Microsoft Corporation.

Software Revision

This guide is valid for A.01.xx revisions of the Agilent G1369A LAN Interface software, where xx refers to minor revisions of the software that do not affect the technical accuracy of this guide.

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 guide contains information to install the LAN Interface (G1369A).

1 Introduction - Around your LAN Interface

In this chapter you will find an introduction to the LAN Interface and its function.

2 Getting Started

In this chapter you will find instructions to help you to set-up your LAN Interface.

3 Getting Help

In this chapter you will find support information about troubleshooting, repair and the Agilent web.

Introduction to the LAN Interface	8
LAN control - what exactly does it do?	9
LAN Interface - what has to be done?	9
LAN control configurations	10
LAN Interface compatibility	12
Installing and cabling the LAN Interface	14
What you will get	14
What you have to do first	15
LAN Interface configuration	18
TCP/IP parameter configuration	18
18	
Configuration switches	19
Initialization mode selection	20
Link configuration selection	23
Automatic configuration with Bootp	24
Configuring the CAG Bootp server program	24
Storing the settings permanently with Bootp program	29
Manual configuration	30
With Telnet	31
With Handheld Controller	34
PC and Agilent ChemStation setup	39
PC Setup for Local Configuration	39
Agilent ChemStation setup	40
Troubleshooting	42
Link Status LEDs	42
Error Messages	43
Repair and Parts Information	44
Agilent Support Information	45
Firmware Update	45

Contents

Reporting of Problems 45

Agilent Web 45

Glossary 46



1

Introduction - Around your LAN Interface

Introduction to the LAN Interface	8
LAN control - what exactly does it do?	9
LAN Interface - what has to be done?	9
LAN control configurations	10
Local configuration using cross-over cable	10
LAN using a HUB and twisted pair cables	10
LAN with existing customer network	11
LAN Interface compatibility	12

In this chapter you will find an introduction to the LAN Interface and its function.



Introduction to the LAN Interface

The LAN Interface (Local Area Network) is the Agilent replacement for the previously used HP JetDirect card in the Agilent 1100 series HPLC modules and the 8453 UV-vis spectrophometer.



Figure 1 LAN Interface

LAN control - what exactly does it do?

In its simplest form...

- control of your instrument and acquires data “remotely” from your desktop (easier access),
- a direct replacement for GP-IB (HP-IB) interface protocol,
- allows your instrument to be placed anywhere on the laboratory/corporate network,
- improves lab “ergonomics” (better organization),

LAN Interface - what has to be done?

- install LAN Interface into the instrument
- install network interface card (NIC) into PC (if not already pre-installed or on-board).
- connect to instrument
 - direct with cross-over cable or
 - to HUB with twisted pair cable
- configure instrument on LAN

LAN control configurations

The basic LAN configurations are shown below.

Local configuration using cross-over cable

The simplest way is a configuration with a single system.

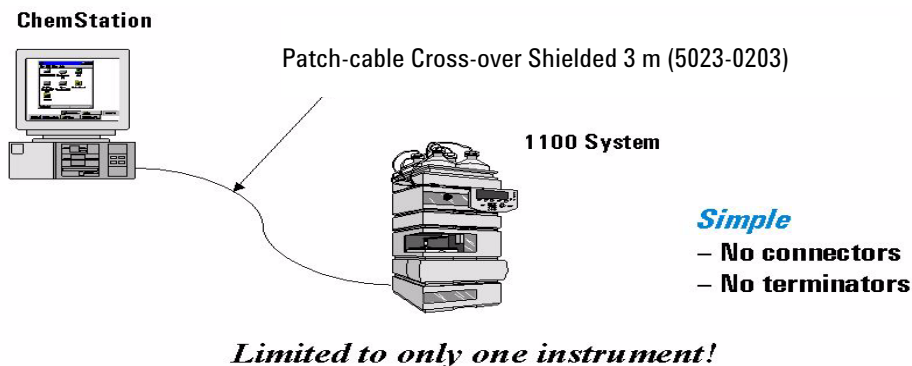


Figure 2 Local configuration using cross-over cable

LAN using a HUB and twisted pair cables

More complicated setup than direct cross-over connection.

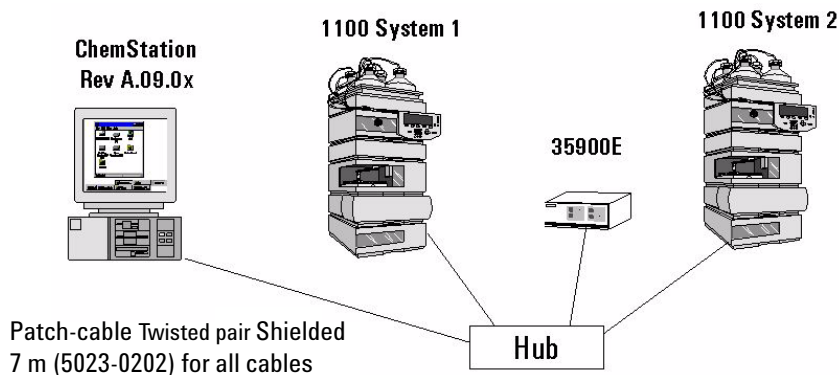


Figure 3 LAN configuration using a HUB and twisted pair cables

LAN with existing customer network

Use MDI/MDI-X port or “Cascade” Port with standard twisted pair cable to connect Hub to a “parent” hub. IP Addresses and other TCP/IP configuration information MUST be provided by the customer’s IT organization. The customer LAN must be able to handle instrument data and must have sufficient bandwidth for instrument acquisition (no overnight backups over the LAN).

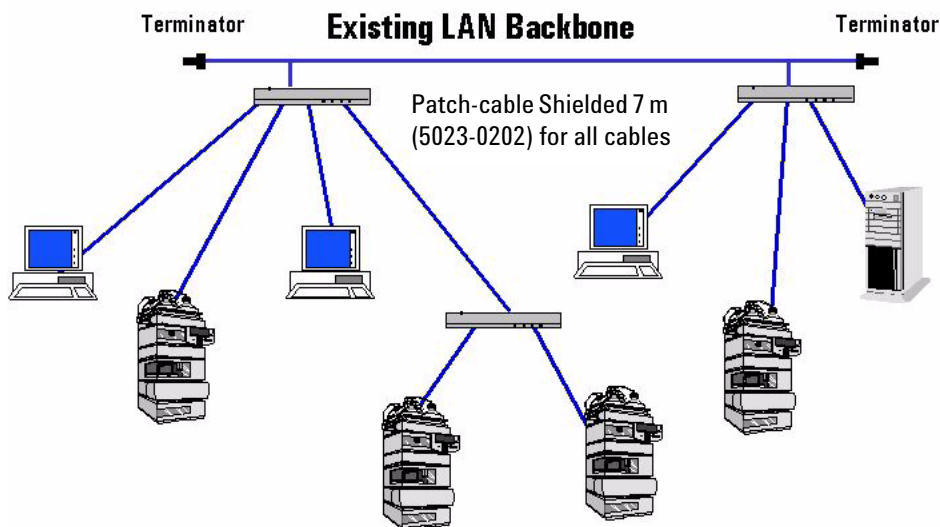


Figure 4 LAN configuration with existing customer network

LAN Interface compatibility

The table below lists the minimum requirements for LAN operation with the LAN Interface.

Table 1 LAN Compatibility

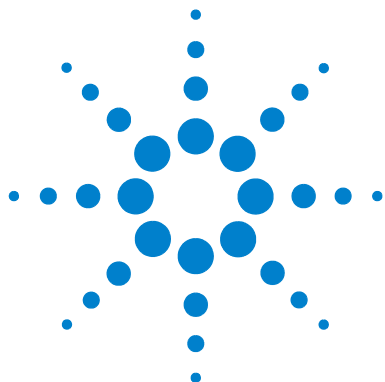
Instrument/Operating Software	Revision (minimum)
Agilent 1100 modules	Firmware A.03.80 and Revision 2 mainboard, see Table 2
Agilent Control Module G1323A	All revisions show just the status page, no editing possible
Agilent Control Module G1323B	All revisions below B.02.02 show just the status page, no editing possible. With B.02.02 and above editing is possible.
Agilent 8453 Spectrophotometer	Firmware 3.30
Agilent Control Module G1818A	No viewing or editing possible
Agilent ChemStation software	A.06.02 or later

LAN compatibility on early 1100 modules

All 1100 Series HPLC modules shipped prior to 1997 are NOT compatible with the LAN Interface communication. The modules which host the LAN Interface (usually the detector module) requires a new main board. The serial number break of the 1100 series modules and the part numbers for the new boards are listed below.

Table 2 LAN compatibility on early 1100 modules

1100 Module	S/N break	P/N Mainboard
G1310A	below DE64300355, US64400233	G1311-66520 or higher
G1311A	below DE64301137, US64401134	G1311-66520 or higher
G1312A	below DE64300703, US64400425	G1312-66520 or higher
G1313A	below DE64302092, US64400886	G1313-66520 or higher
G1314A	below JP64201926	G1314-66521 or higher
G1315A	below DE64301532, US64400333	G1315-66520 or higher



2 Getting Started

Installing and cabling the LAN Interface	14
What you will get	14
What you have to do first	15
LAN Interface configuration	18
TCP/IP parameter configuration	18
Configuration switches	19
Configuration switches	19
Initialization mode selection	20
Using Stored	21
Using Default	22
Bootp	20
Bootp & Store	20
Link configuration selection	23
Automatic configuration with Bootp	24
Configuring the CAG Bootp server program	24
Storing the settings permanently with Bootp program	29
Manual configuration	30
With Telnet	31
With Handheld Controller	34
PC Setup for Local Configuration	39
Agilent ChemStation setup	40

In this chapter you will find instructions to help you to set-up your LAN Interface.

The information is based on the Agilent 1100 series HPLC modules, but can be easily transferred to the 8453 UV-vis spectrophotometer.



Installing and cabling the LAN Interface

What you will get

- G1369A LAN Interface
- LAN cables (for part numbers see [“Repair and Parts Information”](#) on page 44)

Patch-cable
Twisted pair Shielded 7 m
(5023-0202)

CD-ROM with the
manual



Patch-cable
Cross-over Shielded 3 m
(5023-0203)

LAN Interface
(1369-60001)

Figure 5 What you will get (Content of G1369A)

What you have to do first

NOTE

Use an ESD (Electro-Static Discharge) wrist strap when handling electronics. Refer to your instrument manual for details.

- 1 Remove the LAN Interface from it's packaging.

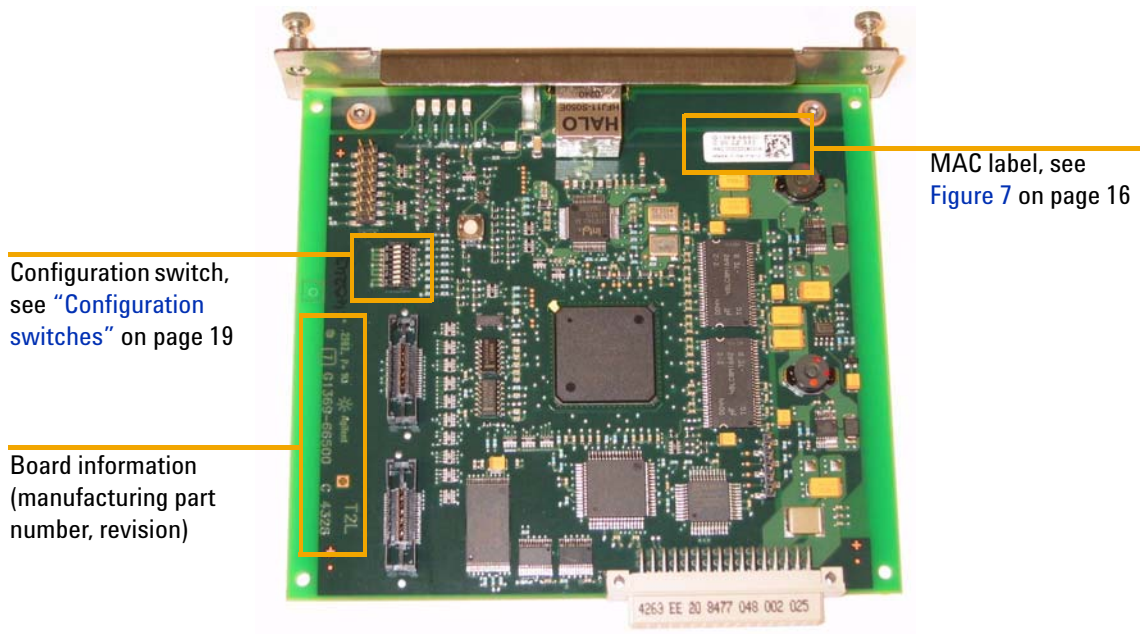


Figure 6 Board Layout

2 Getting Started

Installing and cabling the LAN Interface

- 2 Note the MAC (Media Access Control) address for further reference. The MAC or hardware address of the LAN Interface is a world wide unique identifier. No other network device will have the same hardware address. The MAC address can be found on a label on the card (see [Figure 6](#) on page 15).



Part number of the LAN Interface, see [page 44](#)
Revision Code, Vendor, Year and Week of assembly
MAC address
Country of Origin

Figure 7 MAC-Label

- 3 Turn off instrument line power or remove the power cord before installing the LAN Interface.
- 4 On your instrument, identify the option slot for the MIO accessory card.
- 5 Remove any blank cover plates and ensure that the slot is empty.

NOTE

If the module has the 1100 CAN modification board installed, it probably has a revision 1 mainboard and will not accept the LAN interface. Refer to “[LAN Interface compatibility](#)” on page 12

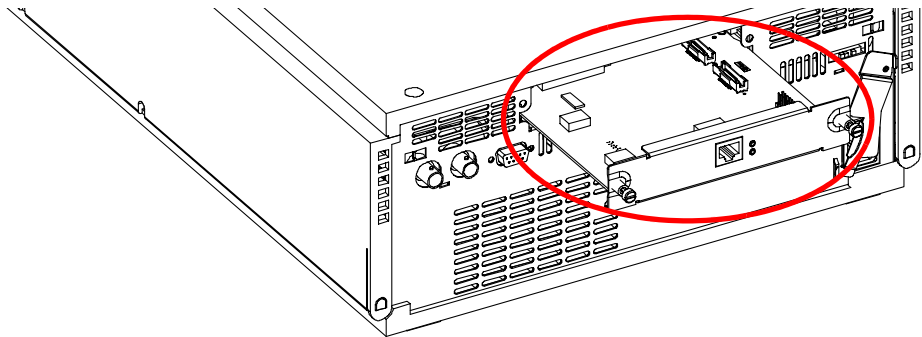


Figure 8 Location of LAN Interface (e.g. 1100 series detector)

NOTE

In 1100 systems, the LAN Interface should be installed in the detector (DAD, MWD, FLD, VWD) due to its higher data handling rate. If no 1100 detector available, use the pump or the autosampler (in this order).

NOTE

The LAN Interface is shipped with the Bootp initialization mode and will use the parameters (IP, Subnet Mask and Default Gateway addresses) from a Bootp server. If you need another initialization mode or other settings, refer to “Initialization mode selection” on page 20 for details before doing the next step.

- 6 Carefully slide the LAN Interface into the slot. Some pressure may be necessary to properly seat the board. Tighten the screws.
- 7 Disconnect your PC from the network and connect the PC network card to the instrument's LAN Interface using a Crossover Network cable (point-to-point) or alternatives, see [page 10](#) and [page 11](#).

CAUTION

Be careful that you connect the LAN cable to the LAN Interface and NOT one of the CAN connections. The CAN bus uses 12-Volt signals, and a misconnection to the CAN bus may destroy network equipment on the other end of the cable.

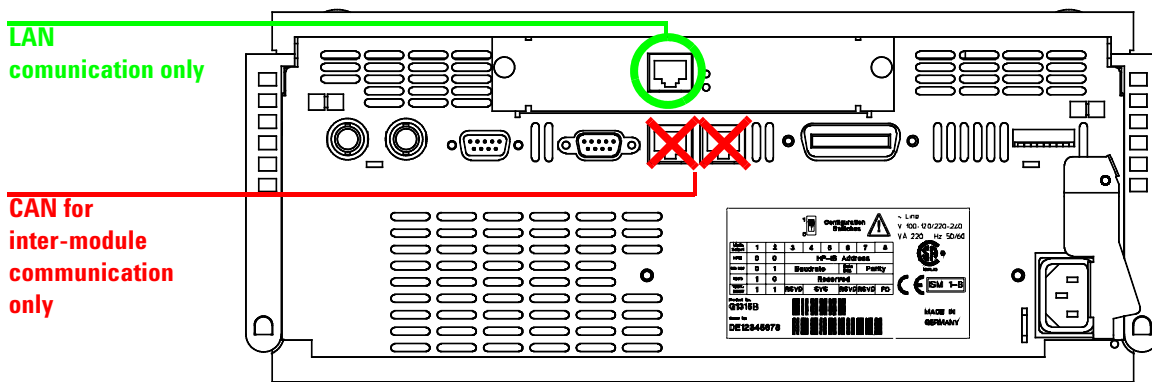


Figure 9 Connect the LAN cable to the correct connector

LAN Interface configuration

TCP/IP parameter configuration

To operate properly in a network environment, the LAN Interface must be configured with valid TCP/IP network parameters. These parameters are:

- IP address
- Subnet Mask
- Default Gateway

The TCP/IP parameters can be configured by the following methods:

- by automatically requesting the parameters from a network-based BOOTP Server (using the so-called Bootstrap Protocol)
- by manually setting the parameters using Telnet
- by manually setting the parameters using the Handheld Controller (G1323A/B)

The LAN Interface differentiates between several initialization modes. The initialization mode (short form 'init mode') defines how to determine the active TCP/IP parameters after power-on. The parameters may be derived from a Bootp cycle, non-volatile memory or initialized with known default values. The initialization mode is selected by the configuration switch, see [Figure 10](#) on page 19.

Configuration switches

The configuration switches are mounted on the card, see [Figure 10](#).

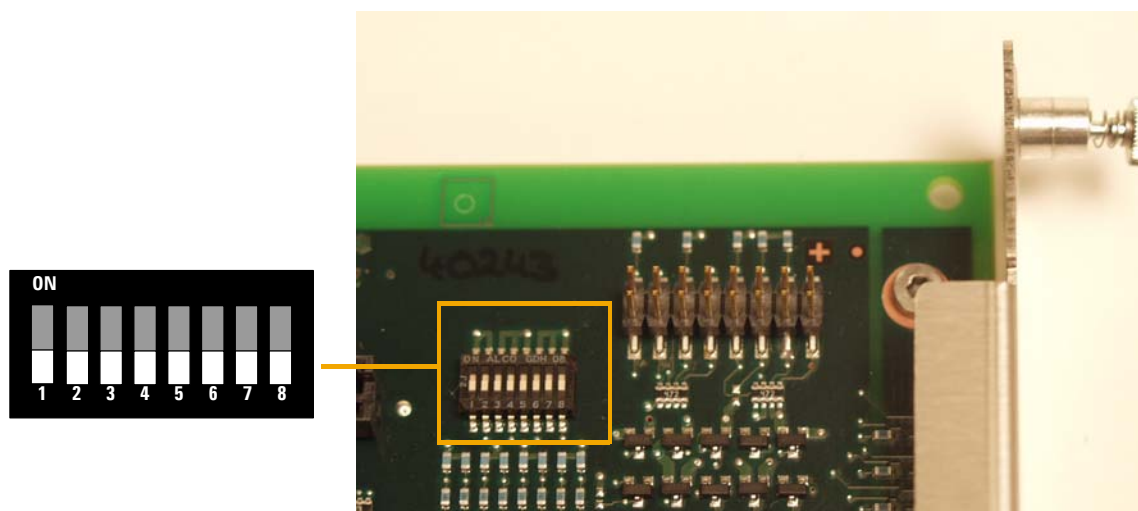


Figure 10 Location of Configuration Switches

The LAN Interface is shipped with all switches set to OFF, as shown above.

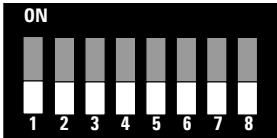
Table 3 Factory Default Settings

Initialization ('Init') Mode	Bootp, for details see page 20
Link Configuration	speed and duplex mode determined by auto-negotiation, for details see page 23

Initialization mode selection

The following initialization (init) modes are selectable:

Table 4 Initialization Mode Switches

	SW 4	SW 5	SW 6	Init Mode
	OFF	OFF	OFF	Bootp
	OFF	OFF	ON	Bootp & Store
	OFF	ON	OFF	Using Stored
	OFF	ON	ON	Using Default

Bootp

When the initialization mode “Bootp” is selected, the card tries to download the parameters from a Bootp Server. The parameters obtained become the active parameters immediately. They are not stored to the non-volatile memory of the card. Therefore, the parameters are lost with the next power cycle of the card.

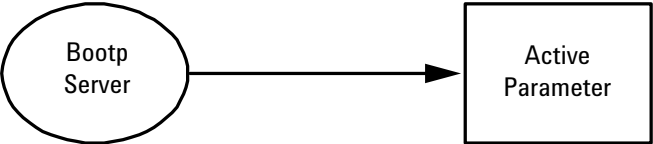


Figure 11 Bootp (Principle)

Bootp & Store

When “Bootp & Store” is selected, the parameters obtained from a Bootp Server become the active parameters immediately. In addition, they are stored to the non-volatile memory of the card. Thus, after a power cycle they are still available. This enables a kind of “bootp once” configuration of the card.

Example: The user may not want to have a Bootp Server be active in his network all the time. But on the other side, he may not have any other configuration method than Bootp. In this case he starts the Bootp Server temporarily, powers on the card using the initialization mode “Bootp &Store”,

waits for the Bootp cycle to be completed, closes the Bootp Server and powers off the card. Then he selects the initialization mode “Using Stored” and powers on the card again. From now on, he is able to establish the TCP/IP connection to the card with the parameters obtained in that single Bootp cycle.

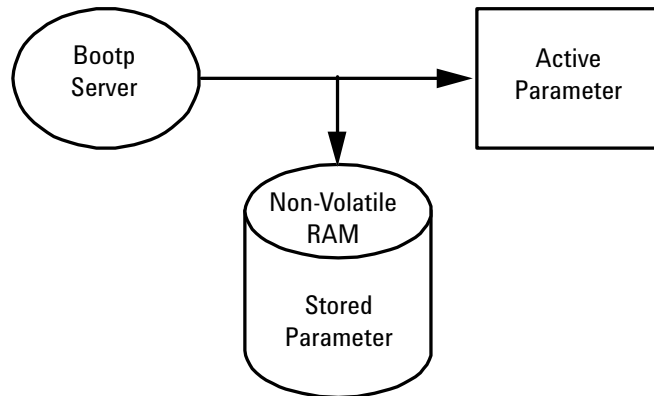


Figure 12 Bootp & Store (Principle)

NOTE

Use the initialization mode “Bootp & Store” carefully, because writing to the non-volatile memory takes time. Therefore, when the card shall obtain its parameters from a Bootp Server every time it is powered on, the recommended initialization mode is “Bootp”!

Using Stored

When initialization mode “Using Stored” is selected, the parameters are taken from the non-volatile memory of the card. The TCP/IP connection will be established using these parameters. The parameters were configured previously by one of the described methods.

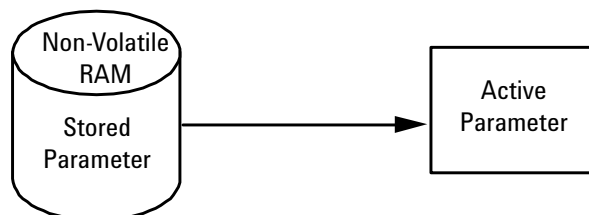


Figure 13 Using Stored (Principle)

Using Default

When “Using Default” is selected, the factory default parameters are taken instead. These parameters enable a TCP/IP connection to the LAN Interface without further configuration, see [Table 5](#).

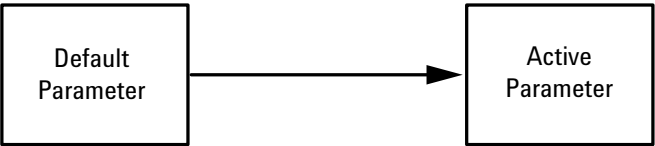


Figure 14 Using Default (Principle)

NOTE

Using the default address in your local area network may result in network problems. Take care and change it to a valid address immediately.

Table 5 Using Default Parameters

IP address:	192.168.254.11
Subnet Mask:	255.255.255.0
Default Gateway	not specified

Since the default IP address is a so-called local address, it will not be routed by any network device. Thus, the PC and the card must reside in the same subnet.

The user may open a Telnet session using the default IP address and change the parameters stored in the non-volatile memory of the card. He may then close the session, select the initialization mode “Using Stored”, power-on again and establish the TCP/IP connection using the new parameters.

When the card is wired to the PC directly (e.g. using a cross-over cable or a local hub), separated from the local area network, the user may simply keep the default parameters to establish the TCP/IP connection.

NOTE

In the “Using Default” mode, the parameters stored in the memory of the card are not cleared automatically. If not changed by the user, they are still available, when switching back to the mode “Using Stored”.

Link configuration selection

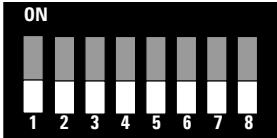
The LAN Interface supports 10 or 100 Mbps operation in full- or half-duplex modes. In most cases, full-duplex is supported when the connecting network device - such as a network switch or hub - supports IEEE 802.3u auto-negotiation specifications.

When connecting to network devices that do not support auto-negotiation, the LAN Interface will configure itself for 10- or 100-Mbps half-duplex operation.

For example, when connected to a non-negotiating 10-Mbps hub, the LAN Interface will be automatically set to operate at 10-Mbps half-duplex.

If the card is not able to connect to the network through auto-negotiation, you can manually set the link operating mode using link configuration switches on the card.

Table 6 Link Configuration Switches

	SW 1	SW 2	SW 3	Link Configuration
	OFF	-	-	speed and duplex mode determined by auto-negotiation
	ON	OFF	OFF	manually set to 10 Mbps, half-duplex
	ON	OFF	ON	manually set to 10 Mbps, full-duplex
	ON	ON	OFF	manually set to 100 Mbps, half-duplex
	ON	ON	ON	manually set to 100 Mbps, full-duplex

Automatic configuration with Bootp

When automatic configuration with Bootp is selected and the LAN Interface is powered on, it broadcasts a BOOTP (Bootstrap Protocol) request that contains its MAC (hardware) address. A BOOTP server daemon searches its database for a matching MAC address, and if successful, sends the corresponding configuration parameters to the card as a BOOTP reply. These parameters become the active TCP/IP parameters immediately and the TCP/IP connection can be established.

Configuring the CAG Bootp server program

NOTE

All examples shown in this chapter will not work in your environment. You need your own IP-, Subnet-Mask- and Gateway addresses.

NOTE

Assure that the LAN Interface configuration switch is set properly. The setting should be either **Bootp** or **Bootp & Store**, see [Table 4](#) on page 20.

NOTE

Assure that the instrument with the LAN Interface installed and connected to the PC is powered off.

NOTE

If the CAG Bootp Server program is not already installed on your PC, then install it from your Agilent ChemStation CD-ROM, located in folder \Bootp.

- 1 The CAG Bootp Server program is placed in the start-up group and automatically is started during the boot process of the PC. It's minimized and located in the task bar.
- 2 Open the Bootp Server window by clicking on it in the task bar.
- 3 Now turn on the module with the LAN Interface and view the Bootp Server window. After some time the Bootp Server will display the request from the LAN Interface with the hardware (MAC) address (this information is also stored in the file trace.txt in the bootp server directory, if Log to Disk is enabled), see [Figure 15](#) on page 25.

The MAC or hardware address of the LAN Interface is a world wide unique identifier. No other network device will have the same hardware address.

The MAC address can be found on a label on the card, see [Figure 6](#) on page 15.

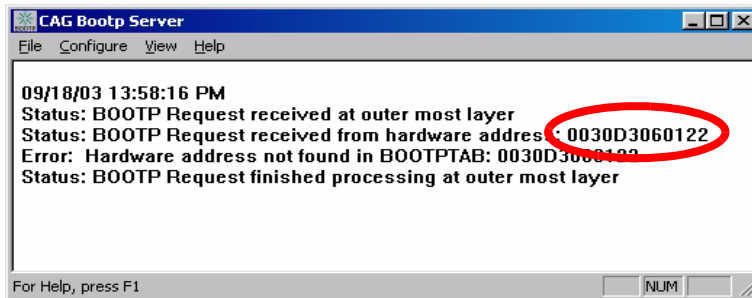


Figure 15 Bootp Server

- 4 Identify your LAN Interface by the MAC address, see [Figure 15](#).

NOTE

If you are working in a network system, you may see other LAN Interfaces appear, overwriting your LAN Interface information periodically.

- 5 Select *Configure* -> *Add Entry* to configure the Bootp Manager ([Figure 17](#)). The drop down box "MAC address" lists all MAC addresses found. Select your MAC address. If no hardware address is found, select *Cancel* and repeat step 3 and step 4.

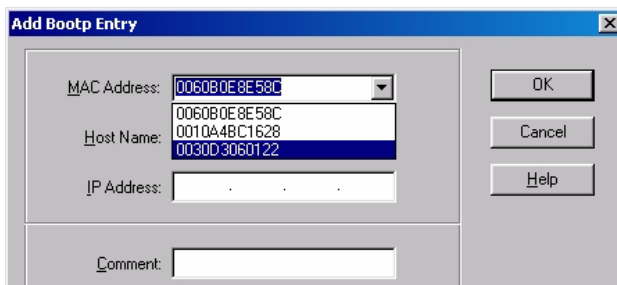


Figure 16 Add Bootp Entry - Select the MAC address

2 Getting Started

Automatic configuration with Bootp

- 6 Specify the Host Name (LC1100-01), the IP address (134.40.24.230), the Comment (LC1100-01) and the Subnet Mask 255.255.248.0 and the Gateway (134.40.24.1).

NOTE

If you are working in a network system, you need your own addresses. Contact your local IT group.

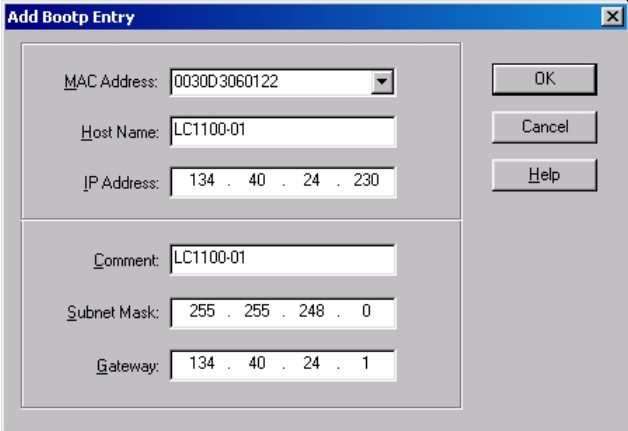


Figure 17 Add Bootp Entry - Enter your parameter

- 7 Exit with *OK*.
- 8 Select *Configure -> Bootp Manager*. All entries made above are shown in [Figure 18](#) on page 27.

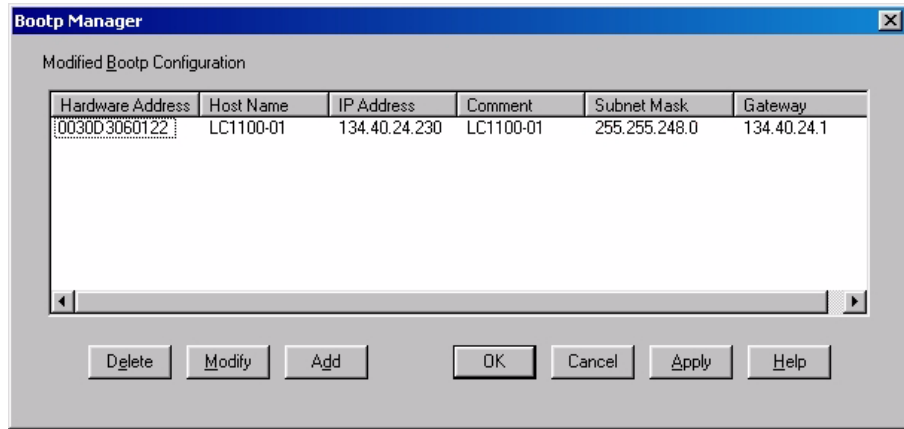


Figure 18 Bootp Manager

- 9 Press *Apply* to activate the changes.
- 10 Press *OK* to exit the Bootp Manager and power cycle the instrument with the LAN Interface, to force it to send a new bootp request again. This time, the MAC address will be recognized by the Bootp Server (Figure 19). It will send the configured IP address and subnet mask information which are necessary for communication to the LAN Interface.

IP address LAN
Interface is
134.40.24.230

IP address PC is
134.40.30.184

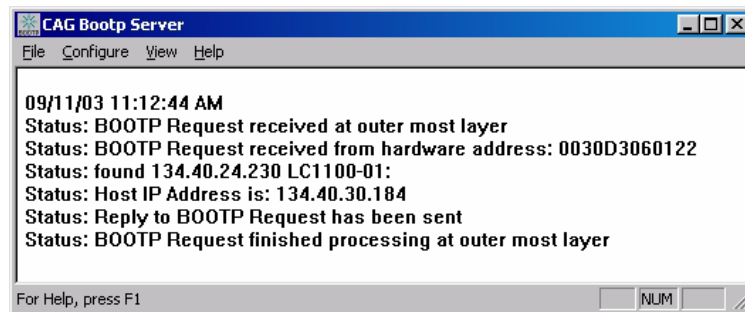


Figure 19 Bootp Server - 1100 module found

2 Getting Started

Automatic configuration with Bootp

NOTE

When using this **Bootp** mode, the parameters are not written into the non-volatile memory of the card. If you delete this Bootp Configuration, the LAN Interface will show up as shown in [Figure 15](#) on page 25 (**Bootp** mode).

If you want to store your parameters permanently on the card (for use without the CAG Bootp server), refer to “[Storing the settings permanently with Bootp program](#)” on page 29.

Storing the settings permanently with Bootp program

If you want to change parameters of the card using the Bootp follow the instructions below.

NOTE

Use an ESD (Electro-Static Discharge) wrist strap when handling electronics. Refer to your instrument manual for details.

- 1 Turn off the module that hosts the LAN Interface and remove the card.
- 2 Change the card's settings of the Configuration Switch to **"Bootp & Store"** mode, see [Table 4](#) on page 20.
- 3 Install the LAN Interface.
- 4 Start the CAG Bootp Server program and open its window.
- 5 If required, modify the parameters for the LAN Interface according to your needs using the existing configuration.
- 6 Press *OK* to exit the Bootp Manager.
- 7 Now turn on the module with the LAN Interface and view the Bootp Server window. After some time the Bootp Server will display the request from the LAN Interface. The parameters are now stored permanently in the non-volatile memory of the card.
- 8 Close the CAG Bootp Server program and turn off the module and remove the LAN Interface.
- 9 Change the settings of the card's Configuration Switch to **"Using Stored"** mode, see [Table 4](#) on page 20.
- 10 Install the card and power cycle the module with the LAN Interface. The card can be accessed now via LAN without the CAG Bootp Server program, refer to ["PC and Agilent ChemStation setup"](#) on page 39.

Manual configuration

Manual configuration only alters the set of parameters stored in the non-volatile memory of the card. It never affects the currently active parameters. Therefore, manual configuration can be done at any time. A power cycle is mandatory to make the stored parameters become the active parameters, given that the initialization mode selection switches are allowing it.

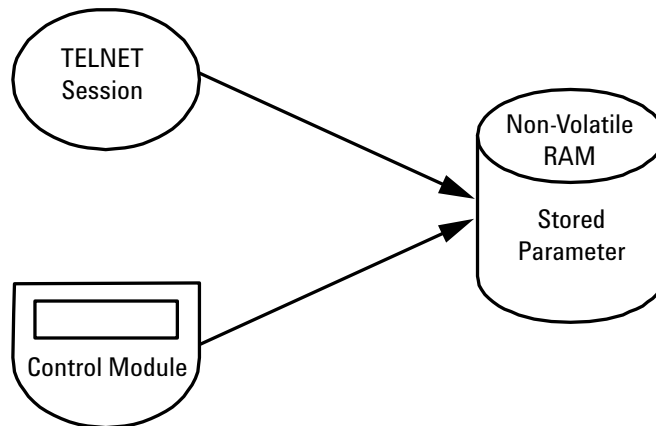


Figure 20 Manual Configuration (Principle)

With Telnet

Whenever a TCP/IP connection to the card is possible (TCP/IP parameters set by any method), the parameters may be altered by opening a Telnet session.

- 1 Open the system (DOS) prompt window by clicking on Windows START button and select “Run...”. Type “cmd” and press OK.
- 2 Type the following at the system (DOS) prompt:

```
c:\>telnet <IP address>
```

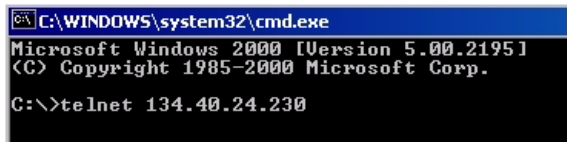


Figure 21 Telnet - Starting a session

where <IP address> may be the assigned address from a Bootp cycle, a configuration session with the Handheld Controller (G1323A/B), or the default IP address (see “[Configuration switches](#)” on page 19).

When the connection was established successfully, the card responds with the following:

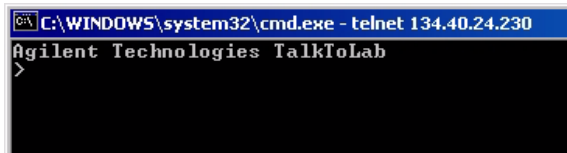


Figure 22 A connection to the module is made

- 3 To change a parameter follows the style:

parameter value

for example: ip 134.40.24.230

then press [Enter], where parameter refers to the configuration parameter you are defining, and value refers to the definitions you are assigning to that parameter. Each parameter entry is followed by a carriage return.

Table 7 Telnet Commands

Value	Description
?	displays syntax and descriptions of commands
/	displays current settings
ip <x.x.x.x>	sets new ip address
sn <x.x.x.x>	set new subnet mask
gw <x.x.x.x>	sets new default gateway
quit	saves changes and exit shell
exit	exits shell without saving changes

NOTE

Any time during the Telnet session you can type “?” then press [Enter] to view available configuration parameters, the correct command format, and a list of additional commands to display.

4 Use the “/” and press Enter to list the current settings.

```
C:\WINDOWS\system32\cmd.exe - telnet 134.40.24.230
Agilent Technologies TalkToLab
>/
Product ID       : G1369A
Firmware Rev.   : A.01.01
MAC Address      : 0030d3060122
-----
Init Mode       : Bootp
Bootp Server    : 134.40.30.184
-----
TCP/IP Properties
- Active -
IP Address      : 134.40.24.230
Subnet Mask     : 255.255.248.0
Default Gateway : 134.40.24.1
- Stored -
IP Address      : 134.40.24.160
Subnet Mask     : 255.255.248.0
Default Gateway : 134.40.24.1
-----
Controller      : not connected
>
```

information about the card
Product id, firmware revision (A.xx.xx are released versions), MAC address, initialization mode

Initialization mode is Bootp
The connected PC/Bootserver is 134.40.24.184

active TCP/IP settings

stored TCP/IP settings in non-volatile memory (not visible if equal to active TCP/IP settings)
connected to PC with controller software (e.g. Agilent ChemStation), here not connected

Figure 23 Telnet - Current settings in Bootp mode

- 5 Change the IP address (in this example 134.40.24.158) and type “/” to list current settings.

```
C:\WINDOWS\system32\cmd.exe - telnet 134.40.24.230
>ip 134.40.24.158
>/
Product ID       : G1369A
Firmware Rev.   : A.01.01
MAC Address      : 0030d3060122
-----
Init Mode        : Bootp
Bootp Server     : 134.40.30.184
-----
TCP/IP Properties
- Active -
IP Address       : 134.40.24.230
Subnet Mask      : 255.255.248.0
Default Gateway  : 134.40.24.1
- Stored -
IP Address       : 134.40.24.160
Subnet Mask      : 255.255.248.0
Default Gateway  : 134.40.24.1
- User -
IP Address       : 134.40.24.158
-----
Controller       : not connected
>_
```

change of TCP/IP setting

Initialization mode is Bootp
The connected PC/Bootserver is 134.40.24.184

active TCP/IP settings

stored TCP/IP settings in non-volatile memory

last user change (not active yet, requires mode “Using Stored” and re-start)

Figure 24 Telnet - Change IP settings

- 6 When you have finished typing the configuration parameters, type:

`quit` and press [Enter] to store the configuration parameters

or

`exit` and press [Enter] to exit without storing parameters.

If the Initialization Mode Switch is changed now to “**Using Stored**” mode, the instrument will take the stored settings when the module is re-booted. In the example above it would be 134.40.24.158 on QUIT and 134.40.24.160 on EXIT.

With Handheld Controller

To configure the TCP/IP parameters before connecting the card to the network, the Handheld Controller (**G1323B with firmware B.02.02 or above for 1100 series modules only**, see LAN Interface compatibility 12.) can be used.

- 1 Press F5 “Views”, select “System” and press the “Enter” key.
- 2 Press F2 “Configure”, select the module where the LAN Interface is installed and press the “Enter” key (Figure 25).

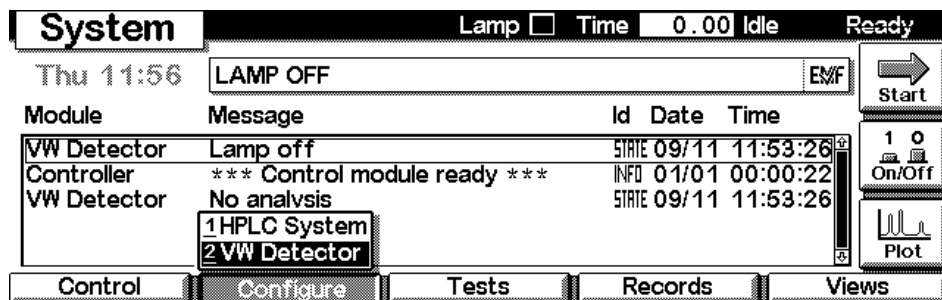


Figure 25 Select module

- 3 Press F1 “Interfaces”, select “MIO” and press the “Enter” key (Figure 26).

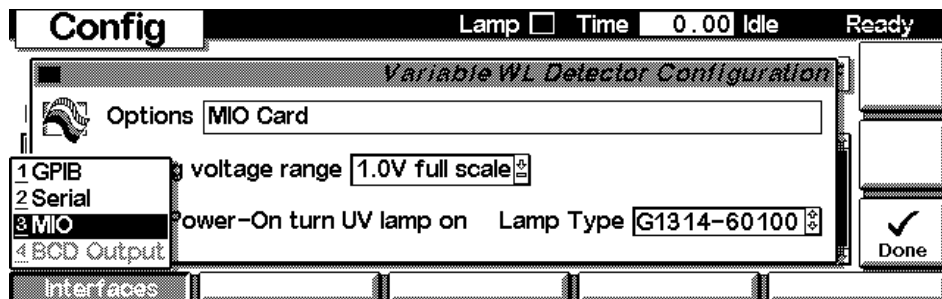


Figure 26 Select MIO

- 4 A Warning message shall pop up. Press “Continue” (Figure 27).

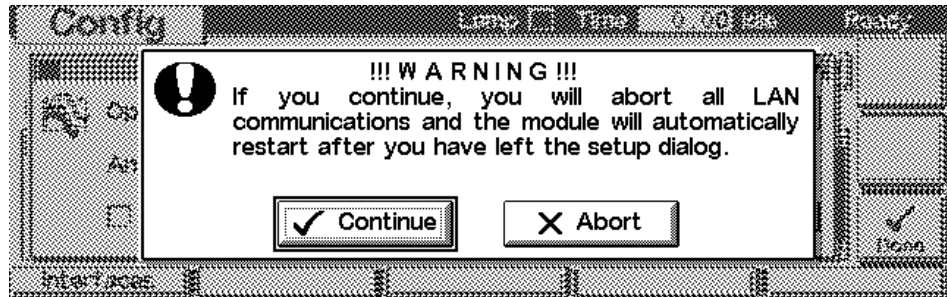


Figure 27 Warning message

- 5 After the Handheld Controller was reading out the LAN Interface you will get an overview of all the parameters that are set in the card (LAN Interface Status Page). The information corresponds to the information in Figure 23 on page 32.

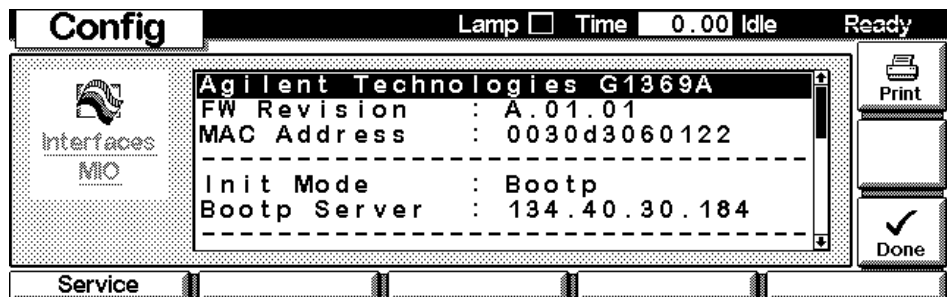


Figure 28 LAN Interface Status Page

In Figure 29 on page 36 the complete listing is shown. For explanations refer to Figure 23 on page 32.

2 Getting Started

Manual configuration

```
Agilent Technologies G1369A
FW Revision   : A.01.01
MAC Address   : 0030d3060122
-----
Init Mode     : Bootp
Bootp Server  : 134.40.30.184
-----
TCP/IP Properties
- active -
IP Address    : 134.40.24.230
Subnet Mask   : 255.255.248.0
Def. Gateway  : 134.40.24.1
- stored -
IP Address    : 134.40.24.160
Subnet Mask   : 255.255.248.0
Def. Gateway  : 134.40.24.1
-----
TCP/IP Status : Ready
-----
Controller    : not connected
```

Figure 29 LAN Interface Status Page (complete)

6 To change the TCP/IP settings, press F1 “Service”.

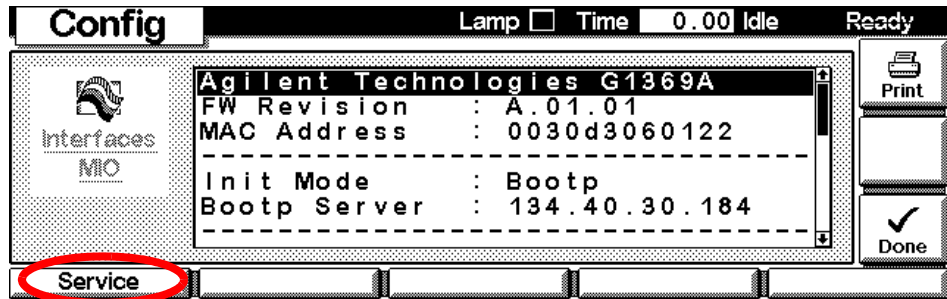


Figure 30 Entering the Service Mode

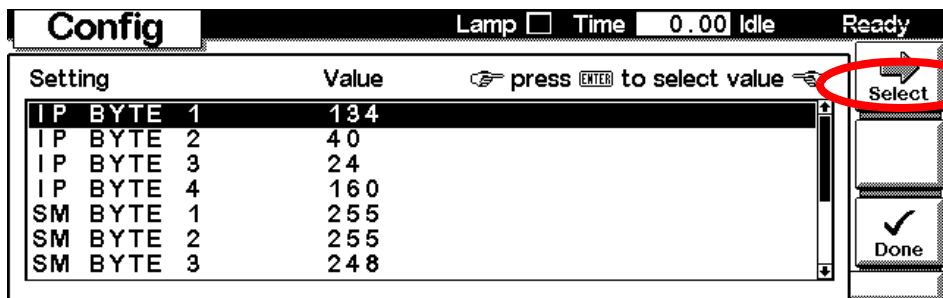


Figure 31 TCP/IP parameters

- 7 Move to the parameter you want to change, enter the new value and press "Enter".

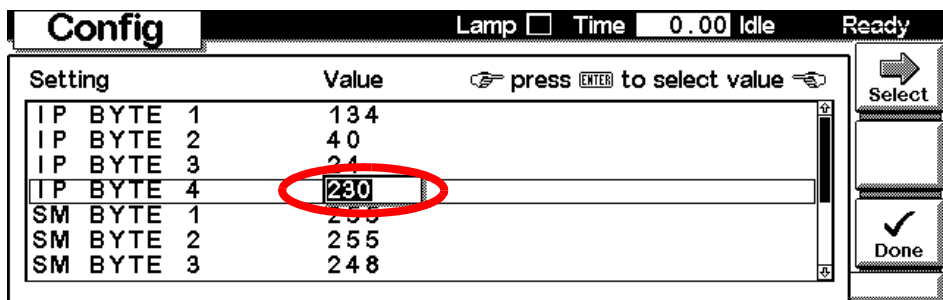


Figure 32 Service

2 Getting Started

Manual configuration

- 8 If you completed your changes, press “Done” to leave the Service section.
- 9 Press F6 “Done” and restart the module by pressing “OK” .

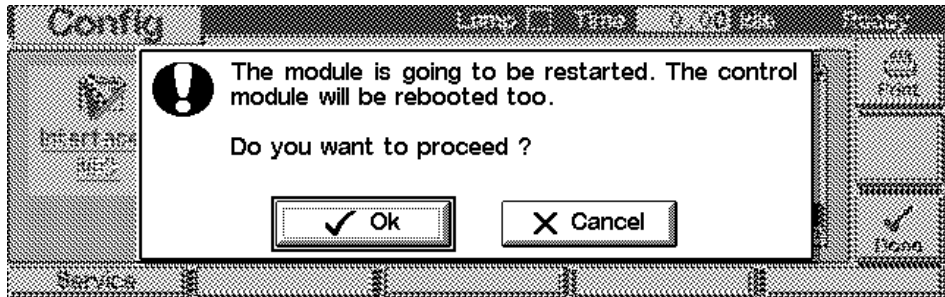


Figure 33 Re-boot screen

PC and Agilent ChemStation setup

PC Setup for Local Configuration

This procedure describes the change of the TCP/IP settings on your PC to match the LAN Interface default parameters in a local configuration (see also “[Local configuration using cross-over cable](#)” on page 10 and “[Using Default](#)” on page 22).

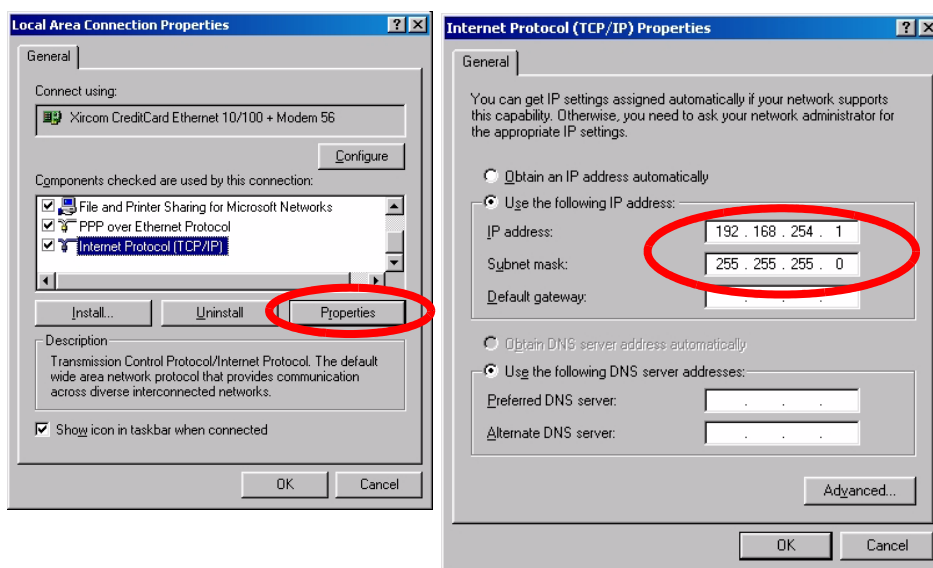


Figure 34 Changing the TCP/IP settings of the PC

Agilent ChemStation setup

- 1 Start the Configuration Editor of the Agilent ChemStation.

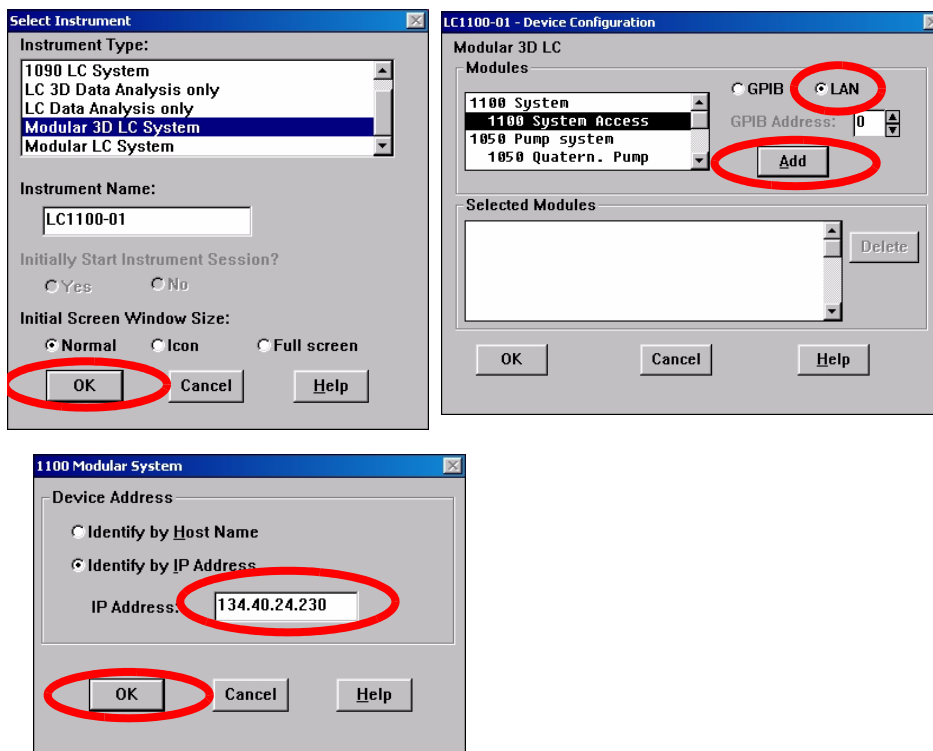


Figure 35 Changing the TCP/IP settings of the Agilent ChemStation

- 2 Add a TCP/IP connection to communicate with the LAN Interface. Use the IP address of the LAN Interface.

NOTE

If using a corporate LAN, IP addresses need to be supplied by the responsible IT department. Also the LAN needs to be able to handle additional traffic.

- 3 Save the configuration, close the Configuration Editor and start the Agilent ChemStation.



3 Getting Help

Troubleshooting	42
Link Status LEDs	42
Error Messages	43
Repair and Parts Information	44
Agilent Support Information	45
Firmware Update	45
Reporting of Problems	45
Agilent Web	45
Glossary	46

In this chapter you will find support information about troubleshooting, repair and the Agilent web.



Troubleshooting

If the LAN Interface does not successfully connect to the network, there are several ways to get status information from the card.

Link Status LEDs

On the card, near the RJ-45 connector, two status LEDs are mounted. See [Figure 36](#).



Figure 36 Status LEDs

The LED named “Speed” shows the actual link speed.

Table 8 LED “Speed”

LED off	link speed 10 Mbps
LED on	link speed 100 Mbps

The LED named “Activity” shows whether the physical link is established or not. In addition, it shows whether the card is transferring data or not.

Table 9 LED “Activity”

LED off	no physical link established
LED on	physical link established
LED blinking	transferring data

Error Messages

The error messages are shown in the LAN Interface Status Page on the Control Module (G1323A/B) only, refer to [“With Handheld Controller”](#) on page 34.

```
Agilent Technologies G1369A
FW Revision   : A.01.01
MAC Address   : 0030d3060122
-----
Init Mode     : Bootp
-----
TCP/IP Properties
- active -
IP Address    : 0.0.0.0
Subnet Mask   : not specified
Def. Gateway  : not specified
- stored -
IP Address    : 134.40.24.160
Subnet Mask   : 255.255.248.0
Def. Gateway  : 134.40.24.1
-----
TCP/IP Status : Error
Bootp timeout
-----
Controller    : not connected
```

Possible reasons:

Bootp server not started or no settings for this MAC address found

Figure 37 LAN Interface Status Page (complete)

If status "Error" shows up, possible error conditions are listed below.

Table 10 Error conditions

Error	Description	Action
Bootp timeout	No reply on Bootp request received	Start Bootp server and/or add settings for the LAN Interface.
Bootp reply incomplete	Bootp reply contained not all information	Complete the necessary information
Gateway in unreachable network	Default Gateway does not match the specified IP address and Subnet Mask	Correct the settings

Repair and Parts Information

The repair level of the product Agilent G1369A LAN Interface is replacement of the complete board.



Figure 38 What you will get (Content of G1369A)

Table 11 Order information

Order number	Description
G1369A	complete product, Agilent G1369A LAN Interface (includes CD-ROM with electronic manual)
G1369-60001	same as G1369A. The board G1369-66500 is a manufacturing number only and cannot be ordered.
G1369-90000	The actual manual as PDF file is available via the Agilent web only, see “ Agilent Web ” on page 45
5023-0203	Cross-over (point-to-point) network cable (shielded, 3 m long)
5023-0202	Twisted pair network cable (shielded, 7 m long)

Agilent Support Information

Firmware Update

The LAN Interface's firmware can be updated, using the firmware provided by the Agilent support web side, see "[Agilent Web](#)". A procedure will be provided with the firmware.

Reporting of Problems

If the LAN Interface shows problems in your system report it with the following information (from the MAC-Label, see [Figure 8](#) on page 16):

- Part number of the LAN Interface
- Board Revision Code, Vendor, Year and Week of assembly
- MAC address
- Installed firmware revision (if known or still accessible, see [Figure 23](#) on page 32 or [Figure 37](#) on page 43).

Agilent Web

Latest documentation or firmware updates for this product (Agilent G1369A LAN Interface) can be obtained from the Agilent web side

<http://www.agilent.com>

> Life Sciences/Chemical Analysis

For firmware select "Technical Support", then look for "Firmware for LC & LC/MS"

For manual select "Library", then search for G1369A and "manual"

Glossary

Table 12 Glossary

Term / Acronym	Definition
10/100Base-TX	Twisted pair Ethernet cable.
Bootp	Bootstrap Protocol, an Internet protocol that enables a diskless workstation to discover its own IP address
CAN	Controller Area Network; a shared broadcast bus, which runs at speeds up to 1Mbit/sec; it is a serial data communications bus for real-time applications.
CAG	Chemical Analysis Group (Agilent term)
DOS	Disk Operating System. The term DOS can refer to any operating system, but it is most often used as a shorthand for MS-DOS (Microsoft disk operating system).
ESD	Electrostatic discharge, the rapid discharge of static electricity from one conductor to another of a different potential. An electrostatic discharge can damage integrated circuits
Ethernet	A local area network (LAN) specified as IEEE 802.3
Gateway	A node on a network that serves as an entrance to another network.
HP-IB or GP-IB	The IEEE-488 Interface Bus (HP-IB) or general purpose interface bus (GP-IB) was developed to provide a means for various instruments and devices to communicate with each other under the direction of one or more master controllers. The HP-IB was originally intended to support a wide range of instruments and devices, from the very fast to the very slow.
IP address	An identifier for a computer or device on a TCP/IP network.
Host	A computer system that is accessed by a user working at a remote location.
Hub	Is some kind of router, which allows clients to connect each other.
LAN	Lab Area Network

Table 12 Glossary

Term / Acronym	Definition
LED	Light Emitting Diode
MAC address	Media Access Control address, a hardware address that uniquely identifies each node of a network.
MIO	Modular Input/Output; interface specification from Hewlett-Packard
RJ-45 connector	Registered Jack-45, an eight-wire connector used commonly to connect computers onto a local-area networks (LAN), especially Ethernets. RJ-45 connectors look similar to the RJ-11 connectors used for connecting telephone equipment, but they are somewhat wider.
Subnet Mask	A mask used to determine what subnet an IP address belongs to. Subnetting enables the network administrator to further divide the host part of the address into two or more subnets.
TCP/IP	Transmission Control Protocol/Internet Protocol; LAN (Ethernet) protocol
Telnet	A terminal emulation program for TCP/IP networks such as the Internet.

In This Book

This guide contains information to install the LAN Interface (G1369A).

- Introduction - Around your LAN Interface
- Getting Started
- Getting Help

© Agilent Technologies, Deutschland GmbH 2003

Printed in Germany
10/2003



G1369-90000



Agilent Technologies

www.agilent.com

© Agilent Technologies, Deutschland GmbH 2003

Printed in Germany
10/2003



G1369-90000



Agilent Technologies

www.agilent.com

© Agilent Technologies, Deutschland GmbH 2003

Printed in Germany
10/2003



G1369-90000



Agilent Technologies