

Agilent 1100 Series Handheld Control Module

User's Guide



Agilent Technologies
Innovating the HP Way

© Copyright Agilent
Technologies 2000

All rights reserved.
Reproduction, adaption,
or translation without
prior written permission
is prohibited, except as
allowed under the
copyright laws.

Part No. G1323-90006

Edition 01/00,

Printed in Germany

Warranty

The information
contained in this
document is subject to
change without notice.

*Agilent Technologies
makes no warranty of
any kind with regard to
this material,
including, but not
limited to, the implied
warranties or
merchantability and
fitness for a particular
purpose.*

Agilent Technologies
shall not be liable for
errors contained herein
or for incidental or
consequential damages
in connection with the
furnishing, performance,
or use of this material.

WARNING

For details of safety,
see Safety Information
on page 142

Warning Symbols Used In This Book



The apparatus is marked
with this symbol when
the user should refer to
the instruction manual
in order to protect the
apparatus against
damage.

User's Guide

In This Book

This book describes how to operate the Agilent 1100 Series modules and systems for HPLC using the control module, shown opposite. The control module provides complete local control and monitoring of a single module or an entire Agilent 1100 Series system. There is no data evaluation in the control module. The control module allows you to do a variety of HPLC tasks including automated sample preparation and injection, isocratic, gradient and multiple method analyses.

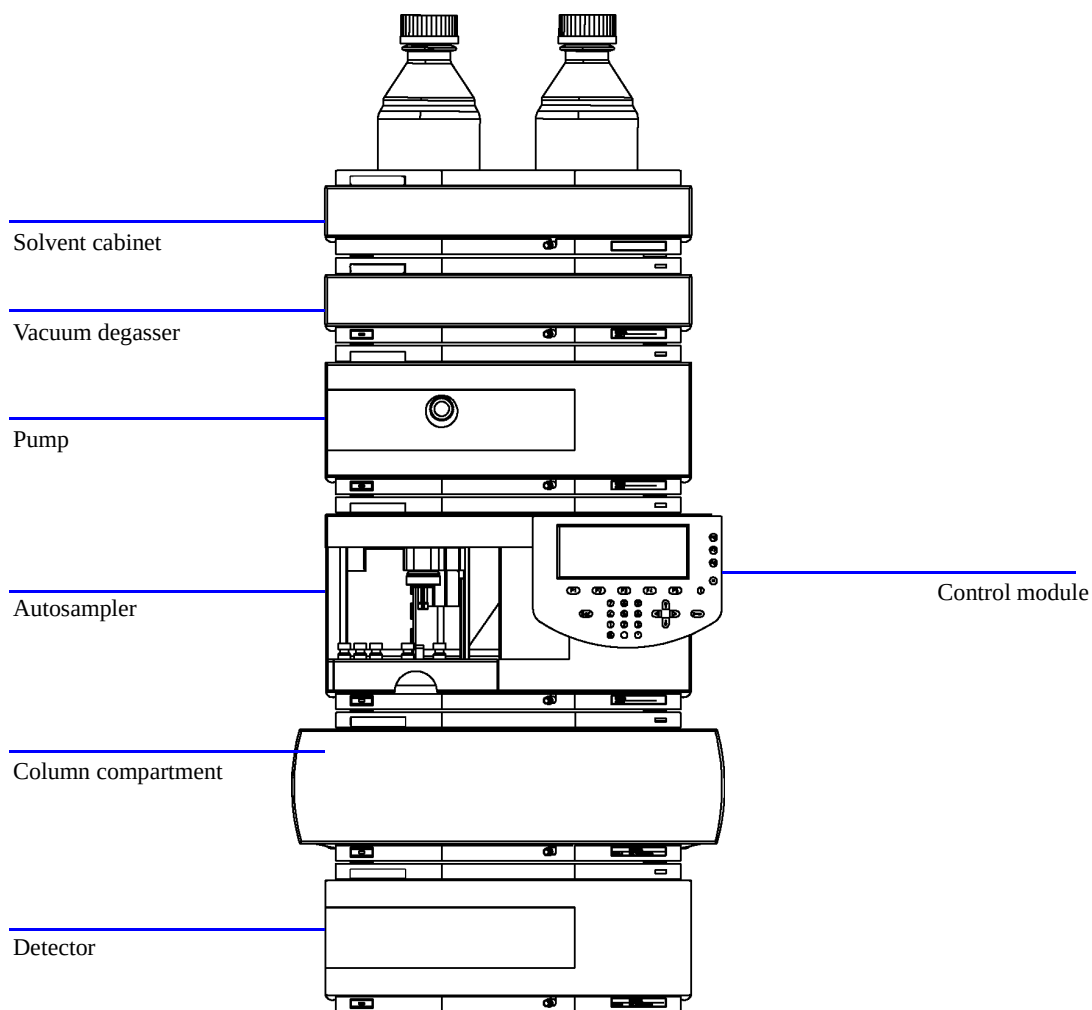
Other Books Available:

Each HPLC module is supplied with a *Reference Manual* which describes installation, troubleshooting, repair, parts and theory of operation.

The control module is supplied with the following:

- *User's Guide*
- *Software Overview Guide*
- *Performance Verification Manual*
- Online Information System

In This Book



In This Book

Part 1 “Using the Agilent 1100 Series Control Module” describes the control module, its features and its functionality.

Part 2 “Using the Agilent 1100 Series Modules” describes how to use the individual HPLC modules to run isocratic, gradient and multiple-vial analyses using a single method or more than one method.

Part 3 “Using the Agilent 1100 Series LC System With Control Module” describes how to run isocratic, gradient and multiple-vial analyses using a single method or more than one method.

In This Book

Part 1 Using the Agilent 1100 Series Control Module

1 The Agilent 1100 Series Control Module

Overview over the Agilent 1100 Series control module 18

Control Module Features 19

Control Module Keys 20

Control Module Software 24

Control Module Versions 31

2 Working with the Control Module

How to use the Agilent 1100 Series control module 33

Installing the Control Module 34

General Functions 35

Troubleshooting 36

Inserting and Removing PC cards 37

Working with Methods 38

Time Programming 43

Automating Analyses 44

Displaying Data Graphically 46

Logbooks 48

Printing Screens 49

Connecting External Devices 51

Firmware 53

Updating Firmware in the Control Module 53

Updating Firmware in the LC Modules 54

Coexecution with Agilent ChemStation 58

Part 2 Using the Agilent 1100 Series Modules

3 Using the Pump

Operational details for the Agilent 1100 Series pumping systems 63

Turning the Pump On and Off 64

Entering Settings 64

Purging the Pump 64

Purging Procedure 65

Adjusting Compressibility 66

Adjusting Stroke Volume 67

Automatic Shut-down 67

Troubleshooting the Pump 68

Tracking Pump History / Pump EMF Limits 68

Resetting the Pump 69

Resetting the Pump Settings 69

4 Using the Degasser

Operational details for the Agilent 1100 Series vacuum degasser 71

Starting the Degasser 72

Removing Gas Bubbles 72

Changing Solvents 73

5 Using the Autosampler

Operational details for the Agilent 1100 Series autosampler 75

Configuring the Autosampler 76

Entering Settings	76
Aligning Transport Arm/Gripper	79
Tracking Autosampler History / EMF limits	79
Resetting the Autosampler	80
Troubleshooting the Autosampler	80

6 Using the Manual Injection Valve

Operational details for the Agilent 1100 Series manual injection valve 81

About the Injection Valve	82
Making An Injection	82
Completely Filling the Sample Loop	83
Partially Filling the Sample Loop	83
How Much Sample Is Actually Injected?	84
How Much Sample Do I Need?	84
An Alternative Way to Fill The Loop	85

7 Using the Detectors

Operational details for the Agilent 1100 Series variable wavelength, multiple wavelength, refractive index, fluorescence light, and diode array detectors 87

Turning the Lamp On and Off	88
Entering Settings	88
Resetting the Baseline	90
Configuring the Detector	90
Troubleshooting the Detector	91
Tracking Detector History	91
Resetting the Detector	92

Lamp Ignition Routine (VWD only) 92
Configuring the Analog Channel Output 92

8 Using the Column Compartment

Operational details for the Agilent 1100 Series thermostatted column compartment 95

Turning the Column Compartment On and Off 96
Entering Settings 96
Configuring the Column Compartment 97
Configuring the Column ID Module 97
Selecting Separated or Combined Mode 98
Selecting the Column Switching Valve (optional) 98
Troubleshooting the Column Compartment 99
Tracking Column Compartment History 99
Tracking Column ID History 99
Resetting the Column Compartment 100

Part 3 Using the Agilent 1100 Series LC System With Control Module

9 Running an Isocratic Analysis

How to analyze the Agilent Technologies isocratic standard sample using a single injection analysis 103

What You Will Need 104
Preparing the LC System 104
Entering Settings 105

Saving Settings in a Method	106
Observing the Chromatogram	107
Starting the Analysis	108

10 Running a Gradient Analysis

How to analyze the Agilent Technologies isocratic standard sample using a gradient analysis 109

What You Will Need	110
Preparing the LC System	111
Entering Settings	111
Saving Settings in a Method	114
Observing the Chromatogram	114
Starting the Analysis	116

11 Running Multiple-Vial Analyses

How to setup multiple vial analyses using the same method and different methods 117

Analyzing Multiple Vials Using the Same Method	118
Analyzing Multiple Vials Using Different Methods	118
Single-Level Calibration Sequences	120
Multiple-Level Calibration Sequences	121
Recalibrating With the Same Group of Standards	121
Recalibrating With Multiple Groups of Standards	123
Synchronizing Analyses with External Devices	126

12 Running an Injector Program

How to create an injector program 129

Creating an Injector Program 130

Entering Injector Program Settings 130

Saving the Method 131

Safety Information 142

Radio Interference 144

Agilent Technologies on Internet 145

Part 1

- 1 The Agilent 1100 Series Control Module 17
- 2 Working with the Control Module 33

Using the Agilent 1100 Series Control Module

The Agilent 1100 Series Control Module

The Agilent 1100 Series Control Module

Overview over the Agilent 1100 Series control module

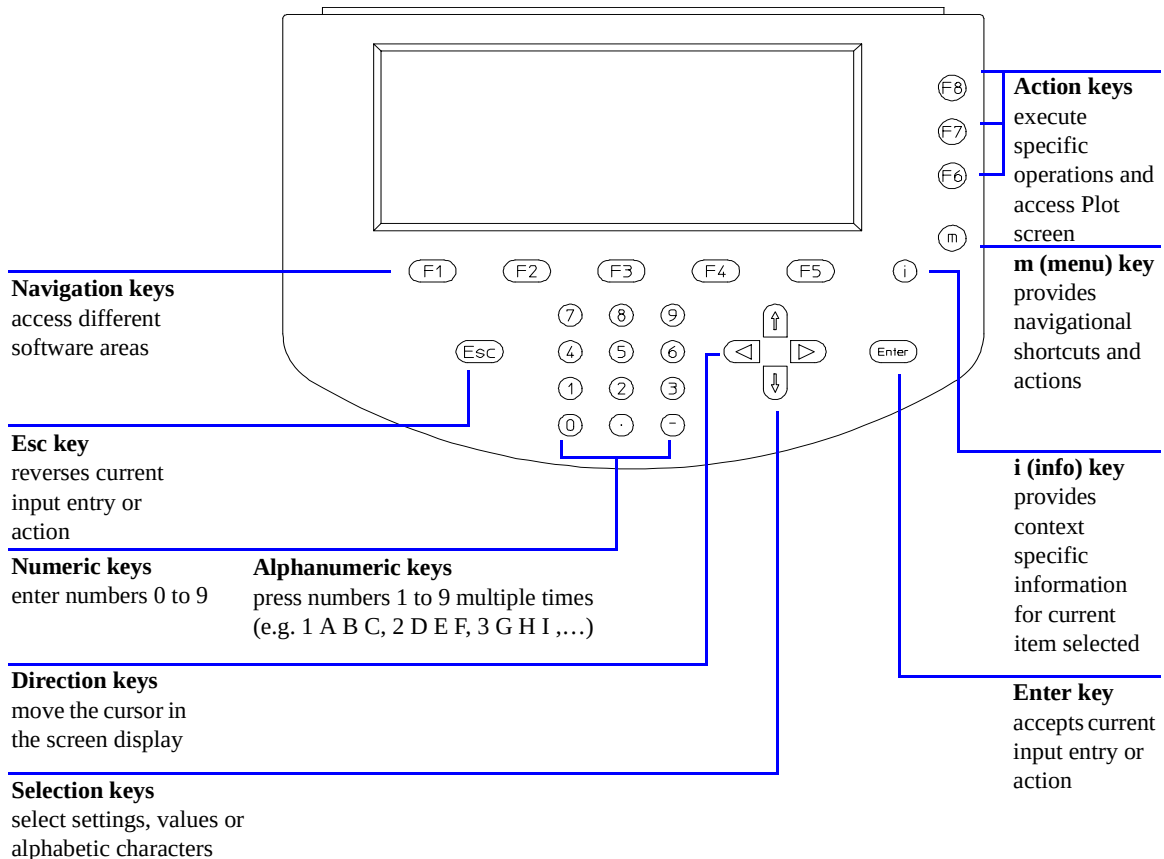
Control Module Features

The control module provides complete local control and monitoring of a single module or an entire Agilent 1100 Series system. You have easy access on every supported function, you can easily control all parameters and settings and you can configure various communication channels with other devices, in order to comfortably analyze the generated data.

- Install any desired configuration of Agilent 1100 Series modules. The control module software will reflect which modules are present in the LC system and adjust the screens accordingly.
- Enter parameter settings for every module, perform reset and on/off functions as well as calibration and configuration settings in a self-explanatory and intuitive way.
- Define automated analyses including methods, timetables, injector programs, method sequences and automated calibration settings using the control module.
- Protect your method from any inadvertent keyboard changes by setting method protection.
- Use PC cards to store and transfer methods and sequences between Agilent1100 systems.
- Monitor all operations and error events using the self-updating logbooks.
- Use the context-sensitive online information system to get further information on all topics.
- Use the context-sensitive menu function to have the quickest access on related functions.
- To help comply with Good Laboratory Practice (GLP) regulations you can select a variety of module tests that will check the performance of the LC system.
- The early maintenance feedback (EMF) limits can be used for scheduling maintenance work.
- Display data graphically using the Plot screen where as many as three different signals can be monitored at the same time.
- Print information to a PCL3 compatible printer connected to the serial RS232 port of a Agilent1100 module.

Control Module Keys

Figure 1 **The Agilent 1100 Series Control Module**



NOTE

A short description of the main keys and the product and serial number is located on the rear of the control module.

Control Module Keys

The display will show you a variety of menu buttons [F1-F5] (in the lower section) or function buttons [F6-F8] (in the right hand section) that can be accessed with the corresponding Navigation (for menus) and Action (for functions) keys.

NOTE

In this context the expression “button” will always refer to a menu or function shown on the display, whereas “key” refers to the actual keys on the keyboard. The key corresponding to a certain button is shown in brackets [F1-F8]

The i (info) key - Online Information System

The online information system provides a quick and convenient way to look up information about a task you are doing or a feature or screen you would like to know more about. The online information system is context-sensitive and provides information related to the current topic.

You can access the online information system by using the i (info) key on the control module keyboard.

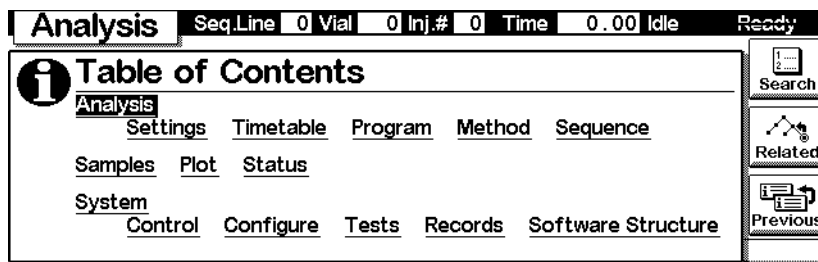
Some words are underlined which indicates available further information. By pressing the Enter key you can display screens providing more specific information about the underlined word.

Some words have a dotted underline which indicates there is an available definition of the word. By pressing the Enter key you can display a small overlay screen with information. You can remove this overlay screen by pressing the Enter key.

- To exit from the online information system press the Esc key.
- To find out what other information topics are available select the Search button [F8]. From there you can choose between Contents and Index.

Figure 2

Online Information System - Table of Contents



- To access further screens containing related information on the currently selected screen select the Related button [F7].

Control Module Keys

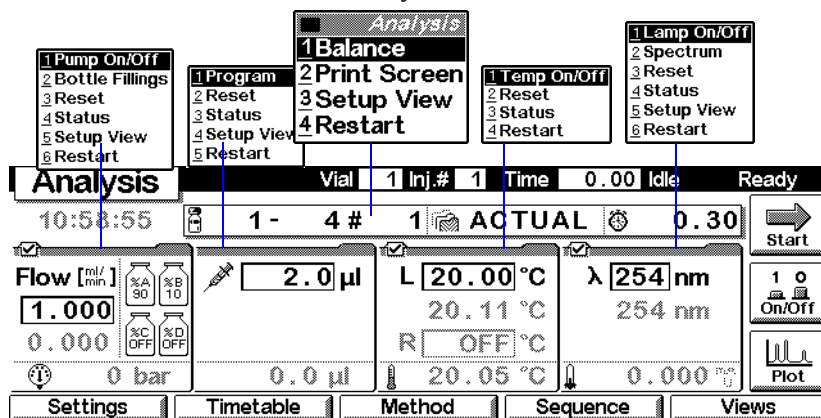
- To move back to the previous information screen select the Previous button [F6].

The m (menu) key

You can use the m (menu) key on the control module keyboard to access selected functionality quickly wherever you are in the software. For example, from the Table of Contents you have quick access to the Print function. In other screens you can easily select the Restart or Default functions or a schematic diagram for explanation purposes. Press the Esc key to exit the menu. See the Quick Reference Guide on the availability of context sensitive menus. In Figure 3 Analysis Screen context menus are shown. Depending on the active parameter entry field context sensitive menus provide different options.

Figure 3

Context Menus in the Analysis Screen



The Esc key

The Esc (Escape) key allows you to exit the current window or screen and leads you back to the last window or screen you were working with.

If you are in one of the main screens, you can use the Esc key to toggle between the current and the previous screen.

In an edit field the previous value can be restored by pressing the ESC key.

The Enter key

With the Enter key you accept a current entry or action. When entering a parameter into a certain field, the Enter key leads you on to the next accessible entry field. In this case it has the same function as the right Direction key.

Navigation keys

These 5 keys [F1 - F5] allow you to switch between the menus. Within these menus the relevant parameters can be set and certain functions can be accessed. The Navigation keys always correspond to a button displayed above them on the screen. The menus accessed via the buttons vary according to the screen you are working with. In some cases pressing a button will cause a list box to appear. From there you must make a choice in order to proceed.

Action keys

The 3 Action keys [F6-F8] trigger a variety of functions. The available functions depend on the screen you are working with.

Selection keys

With the Selection keys (arrow up/down) you can select settings in various list boxes. You can also change values in certain parameter entry fields or enter alphanumeric characters.

Direction keys

With the Direction keys (arrow left/right) you can move back and forth between the entry fields.

Numeric/Alphanumeric keys

These keys allow you to enter numeric values in parameter entry fields. In certain fields where alphabetical characters may be entered you can use the Numeric/Alphanumeric keys to do so. Pressing them several times in sequence changes the current value according to Figure 4.

Figure 4

Keypad of the Control Module

⑦	⑧	⑨
STU	VWX	YZ_
④	⑤	⑥
JKL	MNO	PQR
①	②	③
ABC	DEF	GHI
①	②	③
!/?~	, ; :	+* /

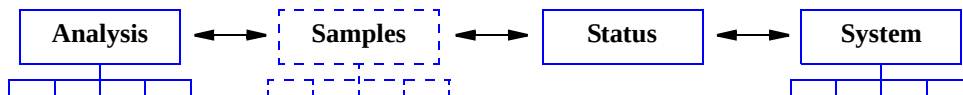
Control Module Software

The Control Module User Interface

In a basic configuration of modules there are at least three main screens, the Analysis, Status and System screens. You can move between these screens using the Views button [F5]. If an Autosampler is part of your system, a fourth screen called Samples can be accessed. Another important screen is the Plot screen which is accessible via an Action button [F6-F8], depending on the active screen.

Figure 5

Main Screens of the Control Module



From these screens other menus can be accessed. When a certain menu is selected by using a Navigation key, in some cases the control module user interface prompts you with a pop-up menu before proceeding. From there you have to make a choice in order to go on, either by using the up/down Selection keys or by pressing the corresponding number on the Numeric/Alphanumeric keys. The latter will give you quicker access to the choices in most cases. Mostly these pop-up menus will

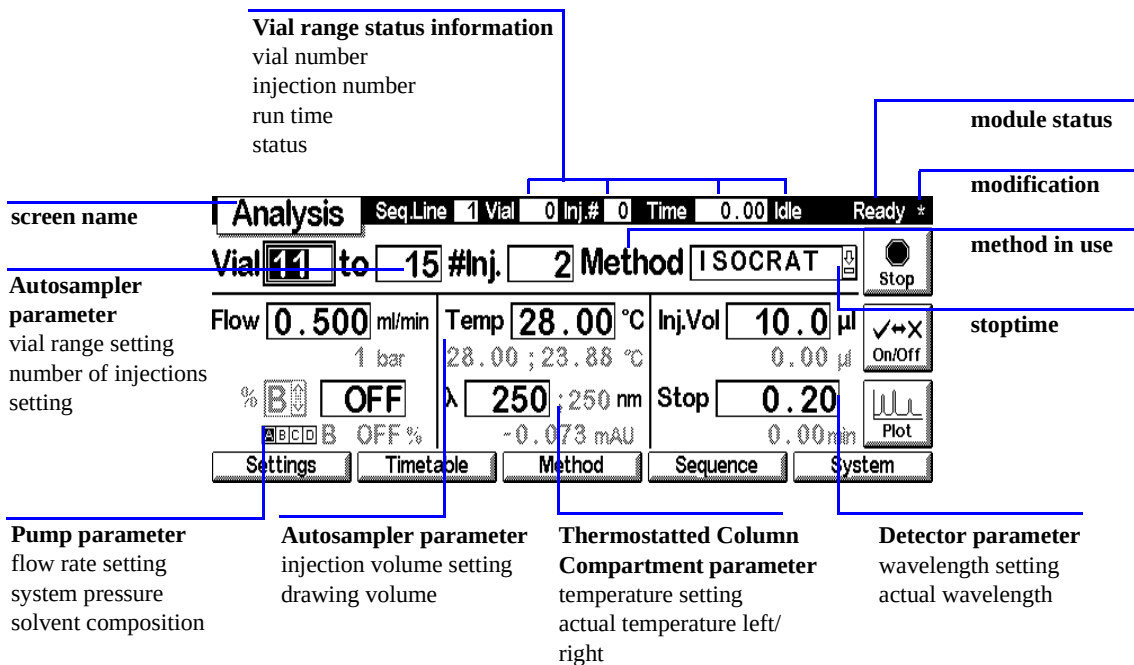
distinguish between the different modules in your system, or in the case of the Views button [F5], between screens.

Analysis Screen

The Analysis screen provides access to all LC analytical settings. The most commonly used settings are displayed with their corresponding set and actual values. Since space on the display is limited you will only see a selection of important parameters. The screen layout will depend on the modules that are included in the system and on the modules selection the user did in the Setup View dialog accessible via the m key (the maximum of visible modules on this screen is 4). The other less used settings of the current method appear in further screens to be accessed from the Settings button [F1]. The actual values are updated continuously

Figure 6

.Analysis Screen Information



The screen provides access to:

- Settings — all LC analytical system and module settings,
- Timetable — time programmable settings,
- Method — method management functions (load, save, delete),
- Sequence — multiple method analyses, and
- Views — other screens: Samples (easy access on vials), Status (most important parameters during analysis) and System (control, configuration, tests and records)

The Start/Stop button [F8] starts or stops a single injection or a multiple injection analysis.

The On/Off button [F7] turns the pump, detector lamp, heater and thermostatted column on or off.

The Plot button [F6] provides direct access to the Plot screen where all important parameters can be displayed graphically. It is possible to display several parameters at the same time.

Using the m (menu) key in the Analysis screen prompts context sensitive menus depending on the active parameter entry field (see Figure 3 on page 22).

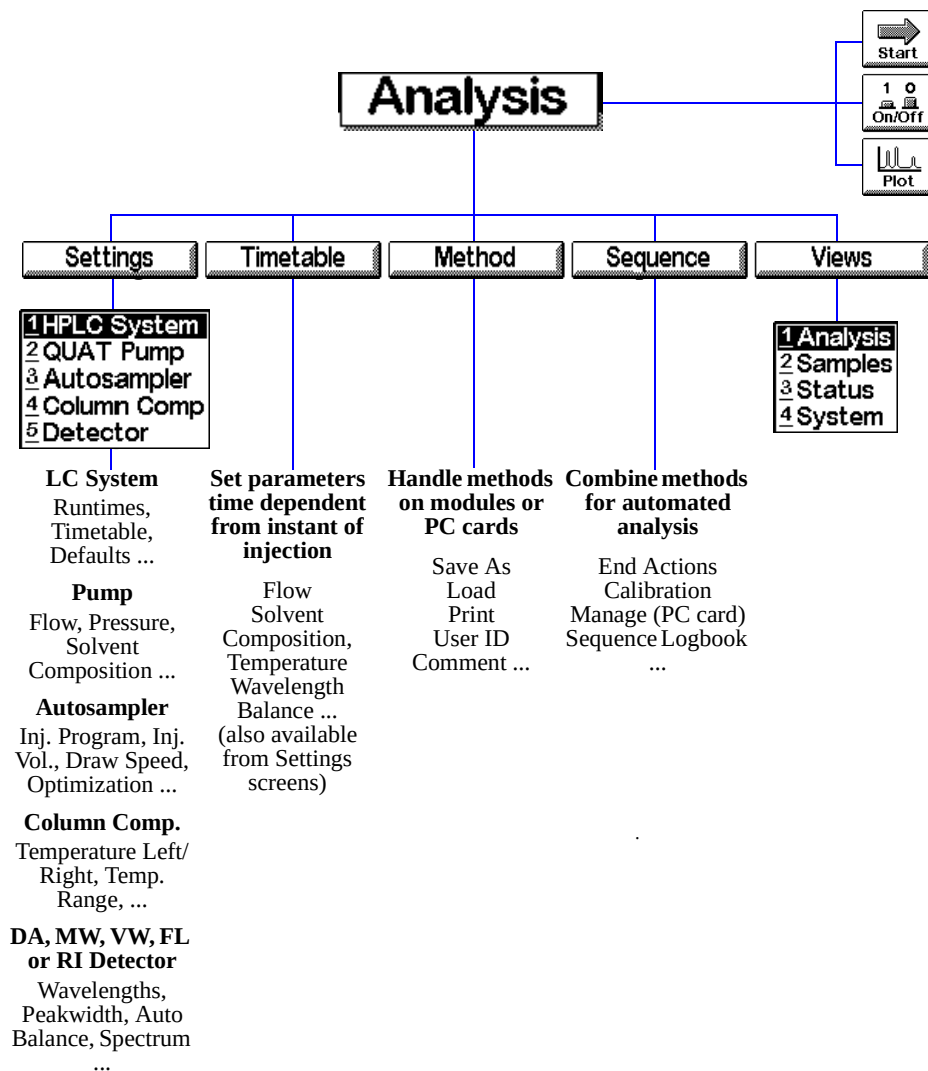
The Module Status on the top right hand side of the screen and on top of the individual module tabs indicates whether the system/ module is ready or not.

Ready indicates the system is ready, **Not Ready** indicates the system/ module is not ready, **ERROR** (only applicable for individual modules) indicates that a vital error has occurred in the module.

The following diagram shows the functionality available in the Analysis screen for an LC system comprising a pump, an autosampler, a thermostatted column compartment and a detector (e.g. Variable Wavelength Detector, Diode Array Detector, Multiple Wavelength Detector).

Figure 7

Where to go from the Analysis Screen - Menu Structure



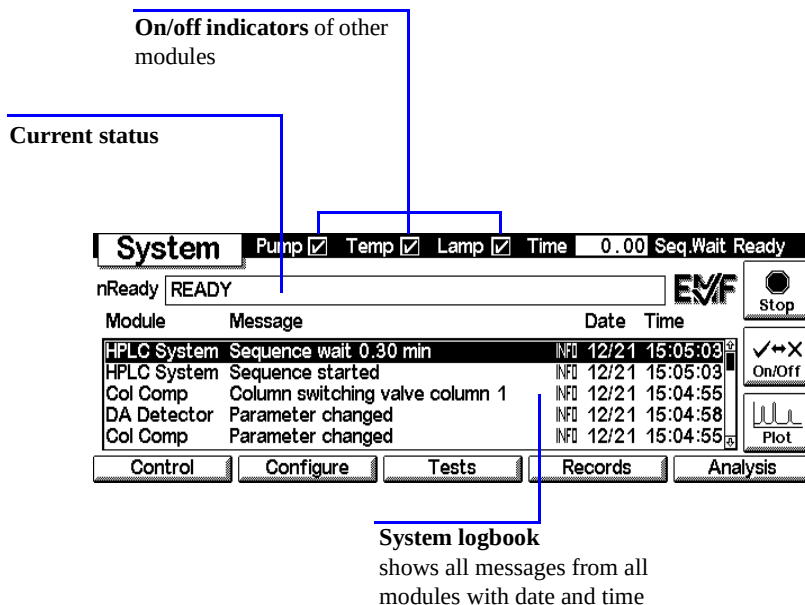
System Screen

The System screen provides access to non-routine settings and provides error checking, control, testing and tracking usage for LC system and modules.

This screen displays the logbook together with the LC system and module status. The EMF (Early Maintenance Feedback) icon will flash to indicate when an EMF limit is exceeded

Figure 8

System Screen Information



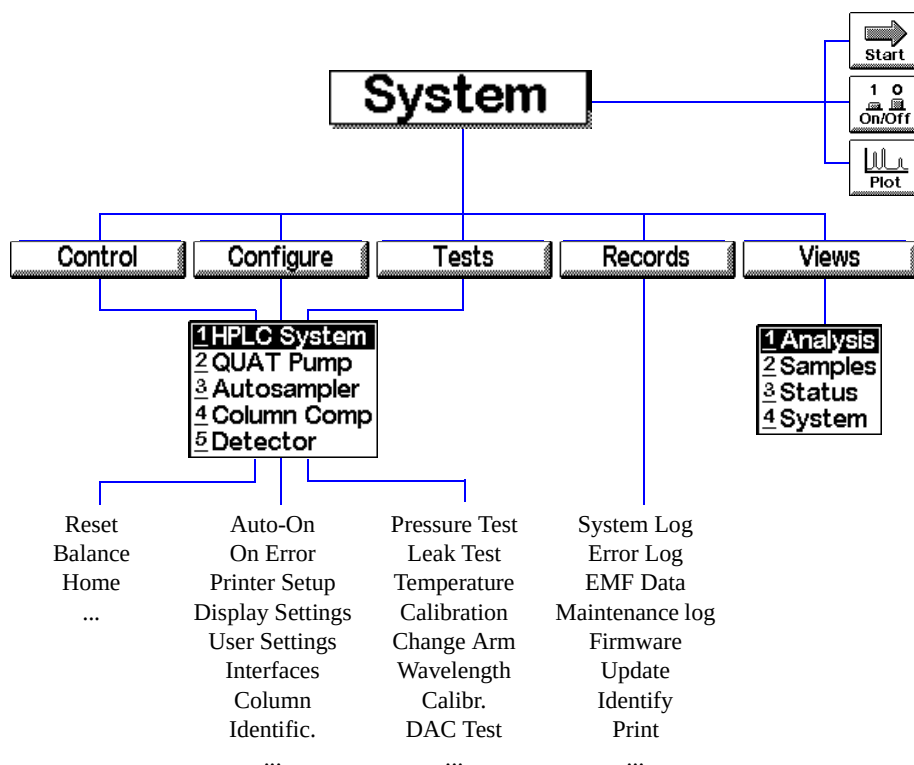
The screen provides access to:

- LC system and module controls, for example, pump on/off, lamp on/off, heater on/off and injector reset,
- LC system and module configuration,
- module tests, and
- LC system and module records.

The following diagram shows the functionality available from the System screen for an LC system comprising a pump, autosampler, diode array detector, variable wavelength detector and thermostatted column compartment.

Figure 9

Where to go from the System Screen - Menu Structure



Status Screen

This screen shows the most important parameters of the system in an easy-to-read style. Included parameters are flow, pressure, column temperature, absorbance units, detector wavelength, elapsed time, etc. depending on the LC system configuration. The screen content is updated in short intervals of roughly one second.

You can also monitor the LC system and module status using the Status bar at the top of each screen. The content of the Status bar varies depending on the screen and is very limited. The status for each module can also be displayed. In the respective Setting screen press the m (menu) key. In the appearing context menu select Status.

Figure 10

Status Screen

Status		Vial	1	Inj.#	1	Time	0.00	Idle	Ready *
Flow	0.000	ml/min	Temp	20.00	°C				
% A	90.0	%		-0.271	mAU				
Pressure	0 bar	Ripple	0.0	%	λ	254	nm		
Elapsed 0.00 min									

Start

1 0
On/Off

Samples Screen

Operating from this screen you have easy access on the sample tray. Using the Vial Range function you can specify the vials you want to analyze and the desired method. A graphical image of the tray offers a good view of the vials chosen. The Sequence option gives you control over the execution of specified analysis sequence steps programmed before. While the analysis is running you can view the processed samples on the tray image.

Figure 11

Samples Screen

Samples		Seq.Line	0	Vial	0	Inj.#	0	Time	0.00	Idle	Ready *
----------------	--	----------	---	------	---	-------	---	------	------	------	---------

☒ **Vial Range** Vial to #Inj

Method Stoptime min

☐ **Sequence** Line to

Start

1 0
On/Off

Plot

End Actions

Timetable

Method

Sequence

Views

NOTE

This screen is only available if an autosampler is part of your system.

Plot Screen

The Plot screen provides an online plot. You can view various signals, depending on the modules in the system, e.g. a chromatogram, pressure signal or temperature signal.

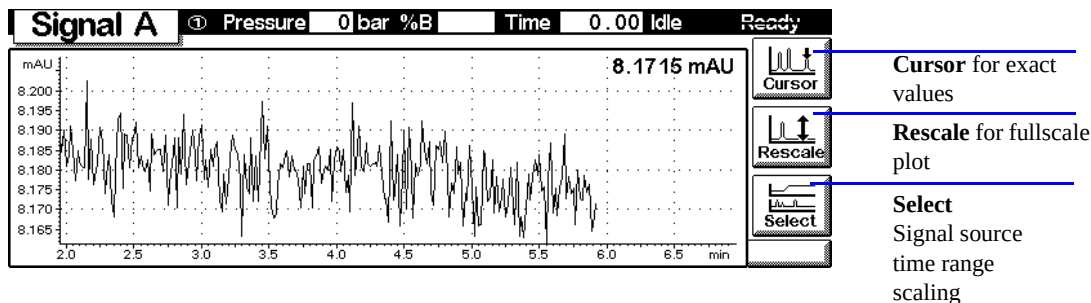
Directly from the Plot screen you can use a cursor to pinpoint information and you can observe information like run time, status, pressure ripple or composition.

Control Module Versions

Using the Direction and Selection keys of the control module you can change the scale of the plot window. With the Rescale button [F7] you can optimize the screen according to the selected signal and parameter ranges.

Figure 12

Plot Screen



Control Module Versions

With the introduction of the Agilent 1100 Multiple Wavelength Detector and the Refractive Index Detector a new Control Module Version, G1323B, was introduced.

To find out which version you possess, check the part number label on the back side of your control module. The G1323A version allows to control the following Agilent1100 modules: Agilent1100 Diode Array Detector, Agilent1100 Fluorescence Detector, Agilent 1100 Variable Wavelength Detector, Agilent 1100 Quaternary, Binary and Isocratic pump, Agilent 1100 Thermostatted Column Compartment, Agilent 1100 Autosampler and Thermostatted Autosampler.

The G1323B version additionally controls the modules Agilent 1100 Refractive Index Detector and Agilent 1100 Multiple Wavelength Detector.

There is no difference in the user interfaces or other functionality.

Working with the Control Module

How to use the Agilent 1100 Series control module

Installing the Control Module

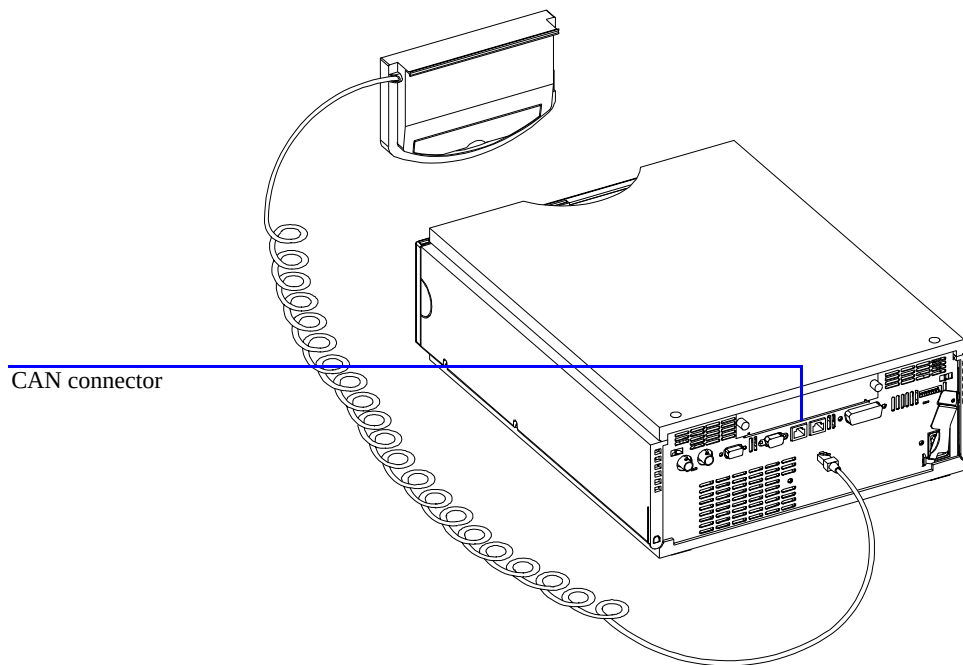
WARNING

The CAN connectors are similar to LAN adapter connectors. Do not insert LAN connectors into the CAN or vice versa, since the CAN uses 24 V and might blow up the LAN card.

Connect the CAN (controller area network) connector of the control module cable to one of the two CAN connectors on one of the Agilent 1100 Series modules.

Figure 13

Connecting the CAN Connector to Rear Panel



General Functions

Turning On/Off LC System and Modules

Use the On/Off button [F7] in the Analysis screen to turn on or off either the complete system or each module individually.

Setting Date and Time

You can set the date and time using the context menu (m key) while the System screen is active. Choose the Date & Time option and press the Enter key. Press the Setup button [F6] to change the settings. Press the Done button [F6] to confirm the new entries. Alternatively, choose Configure and select LC System. Press the Date& Time button [F4]. When date and time are changed in the control module, the settings are automatically stored in all the connected modules. The date and time are backed up by a battery in each module. The control module does not have a clock of its own.

NOTE

Upon startup the modules synchronize their internal clocks. The clocks can also be synchronized by an external chromatographic data system, like the Agilent ChemStation.

Setting Display Contrast

From the System screen choose Configure and select LC System. Press Display and then use the Selection keys to enter values between 0 and 31 to adjust the display contrast according to your personal and location requirements. Then press Done.

Adjusting View

Pressing the m key in the Analysis screen and selecting Setup view allows to select the modules that will be displayed in the Analysis screen. Choosing this option enables you to select the modules present in the Analysis screen. By pressing the Remove/Add buttons [F7, F8] you can move the modules from the Selected Modules to the Available Modules list boxes and vice versa, depending on which module is highlighted.

Configuring the LC System

The LC system is self-configuring to a large extent. It recognizes automatically which modules are installed. The layout of the Analysis screen changes according to the modules present. You can use the Configure button [F2] in the System screen to configure various features such as the Auto-on for the system or Loading (a method), and After Error Condition.

If an error event occurs, you can configure the LC system to load a specified method or turn off the LC system using the After Error Condition setting.

Troubleshooting

Internal diagnostics continuously monitor the module's condition and record any unusual events in an electronic logbook. For example, missing vials or leaking solvent will signal errors and record the errors in the logbook together with the time and date the errors occurred. The logbook is self-updating where the newest entry replaces the oldest entry. If a printer is connected you can easily print out the logbook.

Troubleshooting the Control Module

If your control module does not work correctly, disconnect the module CAN connector from the rear of the Agilent 1100 Series module it is attached to and reconnect it.

If the problem still remains, then

- power down all connected devices and computers and wait 1 minute and then restart, or
- try to use just one Agilent 1100 Series module.

If the problem still remains, call Agilent Technologies.

Inserting and Removing PC cards

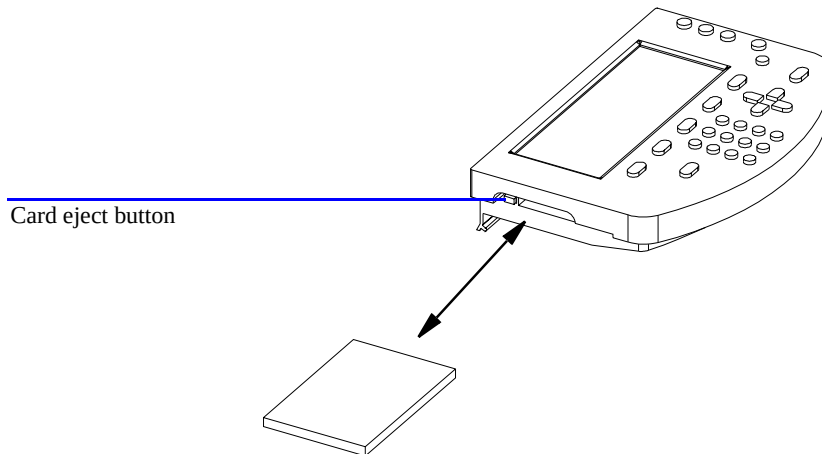
- 1 Insert the connecting side of the PC card into the PC card slot of the control module until you feel it will not go in any further.
- 2 Switch to the Analysis, System or Sample screen, press the m key and select Restart to restart the control module.
- 3 To remove the PC card, press the card eject button on the left hand side of the module.

NOTE

Use only Agilent Technologies or SunDisk Corporation ATA Flashdisks, like the Agilent Technologies F1215C, F1013C or F1014C.

Figure 14

Inserting and Removing the PC card



Working with Methods

A method contains a complete set of injection, separation and detection parameters, including the timetable and injector program. Vial range information is not part of the method.

There are two types of methods:

- The module method. The method parameters are stored in the individual LC modules and not in the control module. A method that is stored in the individual LC modules can be loaded, modified, saved and run from the control module.
- The PC card method. The method parameters are stored on a PC card. A method that is stored on the PC card can be loaded to the LC modules or transferred to another LC system. Methods cannot be run directly from the PC card. The method must first be loaded from the PC card before it can be run. When the PC card method is loaded it becomes a module method.

Unless stated otherwise, the following sections refer to module methods.

Loading a Method

A method can be loaded using the Method button [F3] in the Analysis or Samples screen:

- 1 Enter the Analysis or Samples screen.
- 2 Select Method; the current parameters are displayed.
- 3 Press the Module button [F1].
- 4 Select a method from the list.
- 5 Press the Enter key.

You can also load a method directly from the Analysis screen by using the selection keys and changing the method in the method list.

The Method/Module screen lists all methods that are stored in the modules. For each method there is a date when the method was last changed and a short user description. When a method is loaded it becomes the current method.

A method might be identified as partial method. This means, that there is a mismatch between the actual and the original system configuration, for example

when a module was added to or removed from the system. A partial method can not be loaded as the current method.

Modifying a Method

A method can be modified by changing the settings in the Analysis or Settings screens.

Many of the commonly-used method settings (flow, injection volume, column temperature, wavelength and stoptime) can be modified in the Analysis screen. Other less commonly-used method settings such as eject speed can be modified using the Settings button [F1]. This button displays the Settings menu from which all LC system and module settings can be accessed.

If you change a method setting, the value is immediately downloaded to the LC module. An asterisk (*) will appear in the right corner of the status line to indicate the current method has been modified.

The time-programmable settings can be modified in the Timetable screen.

The injector program settings can be modified in the Injector Program screen accessed from the Autosampler Settings screen.

Specifying a Method Name

- 1 Select the Method button [F3] in the Analysis or Samples screen.
- 2 Select the Save As button [F8].
- 3 Press the Selection key up to enter the letter input mode.
- 4 Select the required letter or number in the method name entry box using the Selection keys. Numbers and the minus sign (-) can be entered directly.

NOTE

Alphanumeric characters can be entered also by pressing the keys 1 to 9 several times (e.g. 1 A B C, 2 D E F, 3 G H I,). See also Figure 4.

- 5 Move the cursor to the next entry position using the right direction key and repeat step 4.
- 6 Repeat step 4 and 5 until you have specified your method name.
- 7 Select the Done button [F6] to accept the method name.

NOTE

The left arrow key can be used as backspace, the right arrow key is used to move to the next character position.

Protecting a Method

To protect the current method.

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select Module from the menu.
- 3 Select the Save As button [F8].
- 4 Select the Protected check box.
- 5 Select the Done button [F6].

The method is now protected against inadvertent changes. Any changes to the method will not be accepted until the method is unprotected, by saving it again without protection. For further protection remove the Control Module from the Instrument and store it in a secure place.

Any unauthorized method or instrument changes can be traced by the system logbook.

Saving a Method

Although it may seem that methods are stored within the control module, in fact all data concerning methods is stored in the modules themselves. The control module generates a list of all available methods that can be loaded.

The number of methods that can be stored depends on the number of timetable and injector program lines included. In general about 15 methods may be stored which contain about 50 timetable lines per module. With differing method contents the actual amount of methods to be stored may change significantly.

Use PC cards in order to store infinite numbers of methods for future use or exchange between LC instruments (see “Transferring Methods Between LC Systems” on page 42).

To save the current method:

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select Save As using the Action keys.
- 3 Enter a name as described in “Specifying a Method Name” on page 39, or simply continue with the next step if you want to keep the current method name.
- 4 Press the Enter key.

NOTE

Optionally you may fill in further data. Activate the Protected check box in order to protect the method. By entering a User ID you can identify your personal modules more quickly. You can also enter a personal comment.

- 5 Select the Done button [F6] to store the method in the modules. Select Yes to confirm save in case you are overwriting an existing method.

The stored method now contains all the current LC system and module settings. The method settings are stored in the individual modules, e.g., all the pump method settings are stored in the pump and not in the control module.

If you disconnect the control module from one LC system and connect it to another LC system, the current method and settings are now shown for the new LC system. To transfer methods from one LC system to another use a PC card.

Deleting a Method

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select Module from the menu.
- 3 Select the method from the method list.
- 4 Select the Delete button [F6].
- 5 Choose Selected Method from the pop up menu. If you want to delete all methods choose All Methods.
- 6 Press the Enter key.
- 7 Select the Yes button to confirm the deletion by pressing the Enter key.

Exchanging Methods With the PC Card

To store and retrieve methods from PC card you can use a comfortable screen that allows easy copying to and from PC card.

Working with Methods

- 1 Insert the PC card into the control module as described in “Inserting and Removing PC cards” on page 37.

NOTE

Methods from a freshly installed PC card are not available. Restart the control module using the Restart function via the m (menu) key from the main screens.

- 2 Select the Method button [F3] in the Analysis or Samples screen.
- 3 Select the PC card button [F2].

NOTE

Use the Initialize button [F1] available in this screen in order to initialize a pre-formatted PC card for use with the control module. This has to take place prior to first use in the control module. Be aware that all data on the PC card will be lost.

- 4 Use the Selection and Direction keys to navigate within and between the method list boxes.
- 5 Use the “Copy >>” and “Copy <<” buttons [F7, F8] to copy methods from and to the PC card.

If you want to store the current method on PC card, you have to use the Save As function (see “Saving a Method” on page 40) in order to store it in the modules first. From there you can copy it to PC card as described in this section.

Transferring Methods Between LC Systems

Methods can be transferred from one LC system to another using a PC card.

- 1 Save the method you want to transfer onto a PC card. See section “Exchanging Methods With the PC Card” on page 41.
- 2 Remove the PC card from the control module.
- 3 Insert the PC card into the control module of the other LC system.

NOTE

If this system does not have a control module connected, use any available control module.

Time Programming

- 4 Restart the control module using the Restart function via the m (menu) key (available from the main screens).
- 5 Load the method from the PC card. See section “Exchanging Methods With the PC Card” on page 41.

Time Programming

To time-program selected settings during the analysis you can create a timetable. Using the Timetable screen, you can create a time based program that will automatically control the pump, detector, column compartment and external contacts.

In some cases the settings will change instantaneously from the initial value to the value specified after a certain time in the timetable (e.g. wavelength). In other cases (solvent composition) these changes take place dynamically, approaching the set value in a step wise and linear manner.

NOTE

The timetable becomes part of the current method when the method is saved.

The timetable is accessed by selecting the Timetable button [F2] in the Analysis screen.

A timetable line can be inserted by pressing the Insert button [F7] and consists of the following:

- Time
Set the time span between the instant of injection and the desired parameter change.
- Module
Choose the module that controls the parameter you want to change.
- Setting
Select the parameter to be changed.
- Value
Enter the desired parameter value.

Automating Analyses

You can edit an existing timetable line by pressing the Enter key. Use the Delete button [F6] and make a choice from the pop-up menu to delete either the selected line or the whole timetable.

You can copy and paste timetable lines by selecting the respective choices available from the context menu (m key).

Automating Analyses

You can use the Sequence screen to create completely automatic unattended analyses from sample preparation to injection. The Sequence screen is accessed by using the Sequence button [F4] in the Analysis or Samples screen.

Using the Sequence screen you can link several methods together. For example, you can first run a method containing an injector program to do sample preparation followed by an analytical run to analyze a batch of samples. You can then run a second method to analyze further samples with different analytical conditions. A delay time can be set in the sequence line. When the second method is loaded, it waits for a specified time before starting the analysis, allowing the column to equilibrate to the new conditions. All sequence events can be traced in the Sequence Logbook available through the Logbook button [F5] in the Sequence screen.

At the end of the sequence you can specify either to load a method (e.g. to flush the LC system to remove buffer salts to avoid crystallization or to program a soft shut-down method) or to turn off the LC system using the End Actions button [F1]. If both options are selected, the shut-down method will be loaded to be available for the next user. However, it will not be executed before the turning off.

You can set up automatic recalibrations using the Calibration Settings screen. This screen allows you to attach calibration settings to a sequence line. The Calibration Settings screen is accessed by selecting the Calibration button in the Sequence screen.

You can recalibrate using one or more standards and have the flexibility of choosing various calibration intervals and patterns. You can define within a sequence line the frequency to recalibrate and the order of calibration vial analysis using the Alter and Multi settings. Alter analyzes the calibration vials alternately. Multi analyzes the calibration vial or vials in complete groups according to the calibration interval.

A sequence line consists of the following:

Automating Analyses

- **Line Number**
Starting with 1 the sequence lines are automatically counted up.
- **Vial Range Information**
Just like in the Analysis screen you can specify a range of vials together with the number of injections per vial.
- **Injection Volume**
Although the injection volume is stored as a method parameter, a sequence has its own injection volume setting (overriding method information). If DEF is specified here the volume as set in the method is kept.
- **Method Name**
Choose a method stored in the modules from the method list box (see “Working with Methods” on page 38). The method has to contain all the relevant parameters as well as timetable or injector program settings.

NOTE

In the Sequence screen you create a program of several methods to be executed in a specified order. Editing of methods or their components (timetables, injector programs etc.) is not possible.

- **Wait Time**
Specify a wait time that creates a gap between method loading and execution. This allows certain module parameters to stabilize before the next analysis is performed.
- **Calibration Settings**
For each sequence line you can define calibration settings by pressing the Calibration button [F1] and choosing Edit/Delete from the pop up menu. You can specify a range of calibration vials, number of injections, injection volume, calibration method and a wait time after method loading. You can also enter specific recalibration parameters, such as the recalibration interval and pattern.

The Online Information System provides accurate information on recalibration options.

Displaying Data Graphically

Using the Plot screen you have many opportunities to display a wide variety of signals on a graphic display while the analysis is performed.

Selecting Signals

Among all the signals available up to 3 can be chosen for graphical display.

- 1 From the Plot screen press the Select button [F6] to show the Plot selection menu.
- 2 Use the Direction and Selection keys to navigate within and between the Available Signals and Selected Signals list boxes.
- 3 Exchange signals between the list boxes by pressing the Move button [F8] or the Enter key.

On the right hand side from the Selected Signals list box you can see the legend to the signals.

You can also enter a time range (X axis) for the plot in this screen.

The different signals can be set up by pressing the Setup button [F7]. Depending on which signal is highlighted you can enter an individual Y-Range setting here.

- 4 When the signals and their X (time) and Y (signal unit) ranges have been specified press the Done button [F6] to switch to the graphic view.

Rescaling the Plot Screen

X (time) axis

To rescale the X (time) axis there are several possibilities:

- Enter a time in the Plot Selection windows (available from the Plot screen via the Select button [F6])
- Perform a rescale directly in the Plot screen by pressing the Direction left/right keys. The right key will shorten the time range by the factor 2. The left key will enlarge the range by the same factor. Press the keys several times to set up the appropriate time frame.

The time range is indicated at the bottom of the Plot screen. This setting is independent from the active signal.

Y (signal unit) axis

To rescale the Y (signal unit) axis there are several possibilities:

- From the Plot Selection windows (available from the Plot screen via the Select button [F6]) choose a signal from either list box and press the Setup button [F7]. You can specify a Y range separately for each signal. This setting can also be made for signals not being part of the Selected Signals list box. Rescaling directly from the Plot screen will overwrite these settings.
- Use the Rescale button [F7] in the Plot screen to adjust the Y axis according to the minimum and the maximum signal value within the set time range. Using this function provides the optimum signal display. It refers only to the active signal indicated at the top of the screen.
- Use the selection keys to change the scaling of the Y axis by a factor of 2 respectively 1/2.

NOTE

Using the m (menu) key and choosing maximize you can enlarge the diagram to full display size. Press the Restore button [F6] to return to the regular view.

Toggling Signals

You can monitor up to 3 different signals from different modules in real-time (updates every second) on the Plot screen. Although all 3 signals are shown in the display, the Rescale button [F7] only refers to the active signal. The active signal is shown in the Plot screen title and can be toggled by pressing the 1, 2, 3 keys on the numeric keypad.

Displaying Exact Signal Values

Selecting the Cursor button [F8] in the Plot screen displays the X and Y value of the current cursor position for the active signal. In this mode, using the Selection keys you can rescale the Y axis. Using the Direction left/right keys you can move the cursor along the graph in an X direction in order to find the position you want.

Print Graph

If a printer is connected to your system you can print the contents of the Plot screen by pressing the m (menu) key and selecting Print Plot. This works also for the maximized view.

Logbooks

The control module keeps track of all kinds of system parameter changes, error messages and maintenance data. You can access the logbooks by pressing the Records button [F4] in the System screen. Information on this screen includes the module product and serial number, firmware version number and the operation time from power-on until now. Additionally an EMF (Early Maintenance Feedback) indicator shows if maintenance is required.

NOTE

You can identify the individual modules by highlighting them in the Records screen and pressing the Identify button [F8]. For several seconds the module's LED will be blinking.

Logbook settings may be entered by pressing the m (menu) key in the System screen and choosing the Setup Logbook option. In the appearing logbook settings menu you can choose which events are to be displayed in the logbooks.

Choose a module from the list box and press the button [F1-F4] corresponding to the type of logbook you want to open.

Maintenance Logbook

On the screen you see the most recent maintenance work together with date and time. Use the Add button [F7] to record maintenance work and update the logbook. The contents of this logbook is permanently stored in the respective modules, but when memory capacity is exceeded the newest entry replaces the oldest entry.

Error Logbook

This logbook displays the most recent error messages with date and time. It is automatically updated whenever an error message is generated by the respective module. This logbook is stored permanently in the module, but when memory is full the newest error will replace the oldest one.

System Logbook

Here all relevant system or event messages are stored until the module is restarted or turned off.

EMF (Early Maintenance Feedback)

In the Records screen the EMF indicator shows if a regular maintenance is needed. The EMF limits can be scheduled by pressing the EMF button [F1] in the Records screen and choosing Setup Limits. Depending on the highlighted module you enter a window where the maintenance intervals can be set. Consult the online information system on Setting EMF Limits.

You can also display EMF events by choosing the Show Events option. This windows shows all EMF limits for the modules that have been operated beyond them. This windows is only updated during startup.

Printing Screens

You can connect a printer with an RS-232 interface to any module using the proper cable.

NOTE

A converter from a serial (RS-232) to a parallel (Centronics) connection is available from Agilent Technologies, part number 5181-1529.

You can configure the printer by pressing the Configure button [F2] in the System screen and then selecting LC System. In the following Config screen press the Printer button [F1] to open the respective dialog. In this dialog you can choose the printer model and the module it is connected to. Define a page layout (including paper size, a banner text and the margins) using the Page button [F8]. Use the Serial button [F7] to enter settings for the RS-232 connection, such as baudrate, bits and parity. This setting can be made individually for each module in the respective configuration settings (see “Connecting External Devices” on page 51).

If the setup is complete you may print a test page using the corresponding button.

NOTE

In many cases a print function is provided via the context menu to be accessed with the m (menu) key, e.g. from the vial range entry fields in the Analysis screen, from the Samples screen and from the online information system.

Print Plot

In the Plot screen you can create a screen hard copy using the m (menu) key and selecting Print Plot from the context menu. The printout will also include a legend and the date and time.

Print Logbooks

Using the m (menu) key to show the context menu in the System screen allows you to choose the Print Logbook function. In the following dialog you can also configure the printer using the Setup button [F8]. The Spooler button [F7] enables you to delete print jobs if several jobs are on the stack. Additionally you may choose the number of copies to be printed.

In the Records screen you can choose the Print button [F6] to create a Configuration Report including the modules' product, serial and firmware version numbers together with the operation time. This button is also available from the System Log, Error Log, Main Log and EMF windows.

Print Method

The current method settings can be printed out using the Print button [F6] in the Method screen. The print dialog as described above will appear. This will print all method settings excluding vial range information.

Print Timetable

Pressing the m (menu) key in the Timetable screen allows you to choose Print Timetable. This will lead to the regular print dialog.

Print Sequence

Entering the Sequence screen and pressing the m (menu) key offers the Print Sequence option.

Print Injector Program

From the Autosampler Settings screen press the Inj. Program button [F3] to enter the Program screen. Via the m (menu) key you have the option to print out the program.

Connecting External Devices

There are several kinds of interfaces that enable the Agilent 1100 Series modules to communicate with a range of other output devices. For some of them extra hardware needs to be installed.

Configuration of selected interface parameters is possible using the Interfaces button [F1] available from the Configure button [F2] in the System screen. This is handled individually for each module, since some interfaces are only available from certain modules (depending on installation).

For further information on interfaces see the corresponding sections in the modules' reference manuals.

APG Remote

Via the 9-pin APG remote connector (included in all modules) the system can communicate with external devices in order to synchronize the analyses. This is necessary when an external device needs some time in order to get ready for a new analysis and thus transmitting of a start request is required (see also "Synchronizing Analyses with External Devices" on page 126 for details on sequence modes). Detailed descriptions of the APG Remote connector are available in the modules' reference manuals.

Among the available signals are:

Power On

This signal is active as soon as all modules connected to the system are switched on.

Shut Down

When the system has a serious problem (e.g. a leak occurs) this alerts all modules to stop relevant operation in order to reduce safety risks.

Stop

This signal asks all modules to reach the ready state as soon as possible. It works only during the analytical run (controlled by the stoptime setting) and causes the system to begin counting down the postrun time.

Connecting External Devices

Ready

When all Agilent 1100 Series modules are ready for the next analysis, this signal is on. Other modules or external devices now can react (e.g. by issuing a start request).

Prepare

This causes the modules to get ready for the next analysis (e.g. the detector will perform a balance).

Start Request

This signal causes the modules to get ready for the analysis (e.g. the autosampler will begin the injection cycle). As soon as all conditions to start the analysis (the injection needle is placed in the seat and the valve is in the proper position) are fulfilled, a Start signal is generated to inform the other modules that now the analytical run starts.

Start

In standard mode only the autosampler creates this signal. This sends an order to start run-time controlled activities to all the modules connected to the APG remote bus. From now on (moment of injection) the runtime counts up.

GPIB

With the GPIB interface (included in all modules) your system is able to communicate with a Personal Computer configured as the Agilent ChemStation. Connect all modules with CAN cables and use one of them to connect to the Agilent ChemStation via an GPIB cable. See “Coexecution with Agilent ChemStation” on page 58 for further details on how to operate the system using the GPIB interface.

Serial / RS-232

Use a standard RS-232 cable to connect a printer to the serial interface (included in all modules). The module communication is enabled with the CAN cables. Choose one module to connect the printer to.

MIO

This interface enables the Agilent 1100 Series modules to communicate with PCs configured as Agilent ChemStations using a local area network (LAN). You can use the MIO interface if the respective extension board is installed in one of your modules and your system is integrated in a LAN.

Firmware

BCD

If the appropriate extension board is part of your system, you can use this output to inform external devices about the vial number currently processed.

External Contacts

With an optional external contacts board you can use various opportunities to synchronize LC activities with external devices.

Firmware

NOTE

When updating firmware using a PC card, make sure the GPIB cable is disconnected from the system and there is only one module (the one where firmware has to be updated) turned on.

The firmware updates can be done using the control module and a PC card. The firmware is loaded from a PC card either into the control module itself or into the modules of the system. You can also update the firmware using a Personal Computer configured as the Agilent ChemStation and connected via the GPIB cable or LAN connection. All Agilent 1100 LC modules can be updated using a Personal Computer and the G1323B control module.

NOTE

The G1323A control module can not be updated using a Personal Computer.

Updating Firmware in the Control Module

- 1 Close the Agilent ChemStation and disconnect the GPIB cable from the system.
- 2 Turn off all modules.
- 3 Insert the PC card into the control module as described in Figure 14 on page 37 .

Working with the Control Module

Updating Firmware in the LC Modules

- 4 Switch on the component where the control module is plugged in.
- 5 In the System screen press Records.
- 6 Choose Controller from the list. If the version number on the display is lower than the current number start the update by pressing FW Update.
- 7 Choose the file with the latest firmware version. For the LC modules file names are 1323A###.BIN where ### is the firmware version number. For the G1323B control module file name are LCB###**.BIN, where ** is the language code, for example EN for English. After selecting the file press Enter. Then press the Execute button [F8] and confirm with Yes to perform the update.

The control module will restart and copy the new firmware into its memory. Firmware loading is indicated by a series of dots displayed on the screen. After reboot the regular (Analysis) screen will be displayed.

In some cases the display contrast has to be readjusted (see “Setting Display Contrast” on page 35).

The correct installation of the firmware can be validated in the System/Records screen. The G1323A/B line should show the new version number.

NOTE

Do not disconnect the cable or power cycle the module while the firmware update is executed.

Updating Firmware in the LC Modules

Within the LC module there are two firmware sections, the resident section (minimum operating system and boot loader) and the operation section. The Transfer button [F7] switches the module from operation firmware to resident firmware mode. In most cases this is the first step in the firmware update process. After the firmware is loaded the transfer command is needed to move the module back from resident to operational mode. Resident firmware is identical for all LC modules whereas operation firmware is different for each module.

NOTE

Update of operation firmware can be done in the resident mode only.

Update of the resident firmware can be done in the normal (operation) mode only.

Updating Firmware of LC Modules via PC

NOTE

The LC modules can also be updated via the firmware upgrade utility delivered with Agilent ChemStation.

Update Functions for Agilent 1100 Series modules

Select File

Allows to select the file on the PC card for the module that is updated with new firmware. Usually binary files are needed:

.BIN for the control module (see “Updating Firmware in the Control Module” on page 53),

.DLB for the module’s operation and resident system.

Transfer

Switches the firmware from the module operation firmware into the resident part. This is the first step in the firmware update process. The resident mode is indicated by a the module’s LED flashing amber. After the firmware is loaded, the transfer command is needed to move the firmware back into the operational part.

Execute

Starts the firmware update. The display shows the progress of the update via the moving bar in the percent field. When updated successfully, “Status: Complete” is displayed.

NOTE

Firmware upgrade of the LC modules resets all EMF counters and erases presently stored methods.

Firmware upgrade of the control module neither resets EMF counters nor affects methods stored in the modules.

Preparing Firmware Update

- 1 Isolate the module you want to update with new firmware by switching off all other modules or by disconnecting all other CAN cables. Remove the GPIB cable from the rear of the module.
- 2 Insert the PC card (as described in “Inserting and Removing PC cards” on page 37) into the control module and reconnect the control module to the selected module to restart the control module.

Switching Firmware Modes

- 1 Prepare the system as described in “Preparing Firmware Update” on page 56.
- 2 Press System — Records and select the module you want to update.
- 3 Press FW Update.
- 4 Press the Transfer button [F7] (and confirm with OK) to switch the selected module into its respective firmware mode.

NOTE

The resident firmware mode is indicated by an amber flashing LED on the module.

After each transfer the control module will reboot. When the resident mode is active the module will be registered as “unknown”. The Analysis screen displays **Resident or Unsupported Module**.

In the System — Records screen you can read the firmware version number of the individual modules. In the normal operation mode you will see the operation firmware version. In resident mode you will see the resident firmware version number.

Updating Resident Firmware

Resident firmware can only be updated in the normal operation mode. Use the Transfer button [F7] as described in “Switching Firmware Modes” on page 56, to check the version number, but be sure to switch back to normal mode afterwards if an update is necessary.

- 1 Prepare the system as described in “Preparing Firmware Update” on page 56.
- 2 Press System — Records and select the module you want to update.
- 3 Press FW Update.

Updating Firmware in the LC Modules

- 4 Select the file for the update (RESXXX.DLB), press the Execute button [F8] and confirm with Yes.

In the screen you can monitor the status of the update procedure.

- 5 When the status bar has reached 100% press the Esc key.

The update of the resident firmware for the respective module has been completed.

Updating Operation Firmware

- 1 Prepare the system as described in “Preparing Firmware Update” on page 56.
- 2 Switch the module you want to update into resident mode as described in “Switching Firmware Modes” on page 56. Also check if a resident firmware update is necessary.
- 3 Press System — Records and select the module you want to update.
- 4 Press the FW Update button [F5].
- 5 Choose the firmware file for the update (####-XXX.DLB), press the Execute button [F8] and confirm with Yes.

NOTE

Make sure you have selected the right firmware and module and the module is in the resident mode as described above.

“Update started” will be displayed on the screen, then the progress of the firmware update can be observed by the moving bar and the % display. If an “update failed, wrong address” error message occurs, make sure that the module is not in the resident mode (indicated by the flashing LED).

- 6 When the operation firmware update is finished press the Esc key to return to the Records screen.
- 7 Press Transfer in the firmware update screen in order to return to normal operation mode (see “Switching Firmware Modes” on page 56).

The control module is rebooted, the status lamp stops flashing and the Analysis screen shows its normal appearance with this module.

- 8 Press System — Records to check that the correct firmware for the module was loaded.
- 9 Reconnect the CAN cables or turn on the other modules.

When updating detector firmware, follow the update procedure with a detector wavelength calibration.

When updating autosampler firmware, make an injection to ensure that sampler alignments are correct. If this is not the case, perform the sampler alignments.

- 10 If other modules' firmware has to be updated, too, repeat the procedure starting with step 1.

Updating Boot Loader Firmware

The Boot Loader firmware (file name BOOT_XXX.DLB) is updated in the same way as resident firmware (see "Updating Resident Firmware" on page 56).

Coexecution with Agilent ChemStation

Features

- Both user interfaces, the control module and the Agilent ChemStation, can be connected to a Agilent 1100 Series system at the same time.
- Parameter entry is possible from both user interfaces. Parameters will be updated on the other user interface within a few moments.
- An Agilent ChemStation sequence can be stopped and aborted from the control module and vice versa.
- The Agilent ChemStation can generate data files from a control module method or sequence. In this case the pre-fix and file name counter in the Single Sample Info section of the Agilent ChemStation must be enabled (protocol mode only).
- If the control module starts an analysis, the Agilent ChemStation is the slave/monitor system.

In general, however, it is not recommended to run both the Agilent ChemStation and the Control Module at the same time. Since the Agilent ChemStation offers a wider variety of controls over the LC Series system and handling it is much more comfortable it should be preferred over the Control Module. Since problems in diagnosis and verification may occur in some configurations the following restrictions apply:

Restrictions

- If a parameter window is open for parameter entry on the Agilent ChemStation,

this specific entry field is disabled on the control module.

- If an analysis is running with the control module, the Agilent ChemStation should not be turned on or rebooted.
- If the Agilent ChemStation starts an analysis, the control module is the slave/monitor system.
- Parameter changes to a method will be identified on the other user interface as modification.
- The control module and the Agilent ChemStation have a different method handling (Agilent ChemStation method can have more information than the method on the control module, e.g. additional DAD parameters that are only accessible from the Agilent ChemStation). To have a method available on both controllers proceed as follows:

If the method is on the Agilent ChemStation and should be saved on the control module or PC card, load the method on Agilent ChemStation and then save the method on the control module (or PC card) with Method — Save As.

If the method is on the control module or PC card and should be saved on the Agilent ChemStation, first load method DEF_LC.M on the Agilent ChemStation (to have no additional parameter in the format) and then load the required method on control module. Then save the method on the Agilent ChemStation with the same name.

WARNING

A method that is available on the control module as protected method can be modified by the Agilent ChemStation and then be saved on the control module without any warning.

Working with the Control Module
Coexecution with Agilent ChemStation

Part 2

3	Using the Pump	63
4	Using the Degasser	71
5	Using the Autosampler	75
6	Using the Manual Injection Valve	81
6	Using the Manual Injection Valve	81
7	Using the Detectors	87
8	Using the Column Compartment	95

Using the Agilent 1100 Series Modules

Using the Pump

Operational details for the Agilent 1100 Series
pumping systems

Turning the Pump On and Off

To turn the pump on or off you can use the On/Off button [F7] in the Analysis or System screen:

- 1 Enter the Analysis or System screen.
- 2 Select the On/Off button [F7].
- 3 Select the Pump button [F8].
- 4 Choose the desired function either with the Action buttons [F6-F8] or by selecting the On/Off buttons in the pop-up window. Uncheck the “to Standby mode only” if you want to turn the pump on or off without going to standby mode.

You can also access this function from the Settings button [F1] available in the Analysis screen or the Control button [F1] in the System screen.

Entering Settings

All pump settings can be entered in the Pump Settings screen. You can access this screen by selecting the Settings button [F1] in the Analysis screen and then choosing the pump. Commonly-used settings such as flow and solvent composition can be set directly in the Analysis screen. The choices available there depend on how many other modules are installed.

Purging the Pump

By opening the purge valve you can redirect the solvent flow from the pump out through the valve exit instead of passing through the injection valve and column. This purging process removes any undegassed or previously used solvent within the pumping system.

Purging Procedure

- 1 Ensure that an outlet tube is connected from the purge valve to a waste solvent bottle.
- 2 Open the purge valve.
- 3 Enter the Analysis screen.
- 4 Set the purge flow rate in the regular flow rate section of the Analysis screen, for example, to 5 ml/min.

NOTE

If a flow rate of more than 5 ml/min is used on a quaternary pump, redefine the upper pressure limit to 200 bar (Accessible from the Analysis screen, press the Settings button [F1], choose pump from the list and press the Enter key. Then choose More ...)

- 5 Set the first channel to be purged to 100%. We recommend you start with the organic solvent channel first. Channel A will automatically be set to 100% when all other channels are set to zero or OFF.
- 6 Turn on the pump using the On/Off button [F7] in the Analysis screen.
- 7 Wait until a continuous stream of solvent comes out of the outlet tube from the purge valve.
- 8 Turn off the pump using the On/Off button [F7] in the Analysis screen and then close the purge valve.
- 9 Repeat the procedure for the other channels you need to purge.

NOTE

The channels of a pumping system are named A, B, C and D (depending on the pump type). %A is automatically calculated by $100\% - (\%B + \%C + \%D)$. If no values for %B, %C and %D are entered, %A is always 100%. To purge the pump you have to go through steps 5 to 8 individually for each of the channels, setting the composition to 100% for the channel to be purged.

Adjusting Compressibility

Liquid chromatographic solvents are compressible under pressure. This solvent compressibility produces a change in flow rate as the pressure changes. To compensate for this effect, you can use the pump compressibility setting to maintain a uniform and accurate flow regardless of the system pressure. The compressibility setting can be accessed by using the Pump Settings button [F1] in the Analysis screen.

Table 1 lists compressibility values for common solvents used in LC.

Table 1

Compressibility Values for Common LC Solvents

Solvent	Compressibility (10^{-6} per bar)
Acetone	126
Acetonitrile	115
Benzene	95
Carbon tetrachloride	110
Chloroform	100
Cyclohexane	118
Ethanol	114
Ethyl acetate	104
Heptane	120
Hexane	150
Isobutanol	100
Isopropanol	100
Methanol	120
1-Propanol	100
Toluene	87
Water	46

When the compressibility setting is set to Off, the pump makes no compensation for the compressibility of the mobile phase.

For each particular compressibility value, the piston stroke (distance piston moves) and the speed at which the piston moves are adjusted accordingly, compensating for the solvent compressibility. For a mixture of solvents we recommend that you choose the compressibility value of the solvent that is present in the highest amount.

Adjusting Stroke Volume

The stroke volume defines the volume of mobile phase which is displaced by one stroke of pump piston 1. You can set the stroke volume to AUTO (automatic) or a value between 20 and 100 μl . When the stroke is set to AUTO, the pump uses large strokes at high flow and shorter strokes at low flow rate. This can improve the mixing performance and gradient linearity by reducing the size of the solvent packets to be mixed.

The stroke volume can be accessed by using the Settings button [F1] in the Analysis screen, selecting the pump and then pressing More ...

The amplitude of the pressure pulsation is directly proportional to the stroke volume. Decreasing the stroke volume decreases the pulsation amplitude, giving a better signal-to-noise ratio with flow-sensitive detectors. To maintain the flow, the pumping frequency is increased, increasing the frequency of the pressure pulsation. Increasing the pump frequency gives better peak area reproducibility at low flow rates.

Automatic Shut-down

At the end of a sequence you can set the pump to be automatically turned off using the Sequence End Actions button [F1] in the Sequence screen. You can turn off the pump completely or load a specific shut-down method.

Troubleshooting the Pump

- 1 Select the Tests button [F3] in the System screen.
- 2 Select Pump from the menu.

You can now select various tests to check the pump. For further information about these tests see the *Reference Manual* for the Agilent 1100 Series pump.

Tracking Pump History / Pump EMF Limits

- 1 Select the Records button [F4] in the System screen.
- 2 Select Pump from the list.
- 3 Press the Enter key.

You can now check the amount of solvent the pump has delivered (liquimeter) and the wear count.

The liquimeter displays the total volume of solvent that the pump has delivered since it was last reset. You can use the liquimeter limit to set up a preventive maintenance schedule for the pump. For example, make a note of the total number of liters pumped when you change the piston seals. This volume is a benchmark, reflecting the lifetime of the seals for your solvents and application. When the limit is exceeded an early maintenance feedback message will appear indicating it is time to change the seals. Changing the seals in advance will prevent you having to repeat analyses, due to leaking seals.

The wear count displays a calculated value indicating the wear of the pump seals. You can also use the seal wear limit to set up a preventive maintenance schedule for changing the pump seals when the limit is exceeded.

Resetting the Pump

- 1 Select the Control button [F1] in the System screen.
 - 2 Select Pump from the menu.
 - 3 Press the Enter key.
 - 4 Select the Reset button [F7] to stop the pump operation and perform a hardware initialization.
-

Resetting the Pump Settings

- 1 Select the Settings button [F1] in the Analysis screen.
- 2 Select Pump from the menu.
- 3 Press the Enter key.
- 4 Select the Default button [F7] to reset the pump settings to their default values.

Using the Pump

Resetting the Pump Settings

Using the Degasser

Operational details for the Agilent 1100 Series
vacuum degasser

Starting the Degasser

- 1 Press the line-power switch on the front panel.
- 2 Turn on your pump by selecting the On/Off button [F8] in the Analysis screen.
- 3 Observe the degasser status lamp on the front panel.

The status lamp is off when there is sufficient vacuum in the degasser.

During operation, the status lamp may be yellow for several seconds. This indicates there is insufficient vacuum in the online degasser and the vacuum pump is on to create sufficient vacuum.

Removing Gas Bubbles

If you see gas bubbles in the solvent tubing or inside the solvent filter:

- 1 Disconnect the solvent tube of the first solvent channel from your pump.
- 2 Connect the syringe adapter onto the syringe from degasser accessory kit.
- 3 Pull the syringe plunger to draw solvent through the degasser and tubing. Continue to draw solvent through tubing until no gas bubbles are visible.
- 4 Disconnect the syringe adapter from the solvent tube.
- 5 Connect the solvent tube to your pump.

Repeat step 1 through step 5 for other solvent channels.

Changing Solvents

If you are changing to a solvent that is immiscible with the solvent currently in the tubing:

- 1 Replace the current solvent:
 - with iso-propanol, if current solvent is organic, or
 - with water, if current solvent is an organic buffer or salt.
- 2 Flush the online degasser and all tubing thoroughly.
- 3 Replace the iso-propanol or water with the new solvent.

Using the Degasser

Changing Solvents

Using the Autosampler

Operational details for the Agilent 1100 Series
autosampler

Configuring the Autosampler

- 1 Select the Configure button [F2] in the System screen.
- 2 Select Autosampler from the menu.
- 3 Press the Enter key.

You can configure the syringe volume, seat capillary volume, action on missing vial, trays, interfaces etc.

Configuring Interfaces

- 1 Select the Configure button [F2] in the System screen.
- 2 Select Autosampler from the menu.
- 3 Press the Enter key.
- 4 Select the Interfaces button [F1].
- 5 Select the interfaces you want to operate (see “Connecting External Devices” on page 51).
- 6 Select the Done button [F6].

Entering Settings

All autosampler settings can be entered in the Autosampler Settings screen. You can access this screen by selecting the Settings button [F1] in the Analysis screen. From here you can access the timetable function (to set up a timetable for the Autosampler only), the runtimes screen and the injector program screen. Commonly-used settings such as injection volume and vial number can be set directly in the Analysis screen.

Entering Settings

More specific settings (draw speed, eject speed, draw position offset) can be entered using the More... button [F1] and choosing the Settings option.

Setting the Injection Mode

There are two modes which can be set:

- Standard

without pre-defined needle wash,

with pre-defined needle wash (wash vial number can be set).

If the pre-defined needle wash function is enabled, the needle moves prior to the injection (with reagent in the needle) into the wash vial to remove sample deposits from the needle outside. This prevents sample stay in the seat and create memory effects on the next injection cycle.

- Injector Program

See “Running an Injector Program” on page 129.

Optimizing Autosampler Performance

There are two optimization modes available on the autosampler (called prefetch and overlap). Both options allow to shorten the analysis time for operations that require a high sample throughput.

The optimization modes can be chosen from the Analysis screen:

- 1 Press the Settings button [F1].
- 2 Choose the Autosampler from the list.
- 3 Select the More... button [F1].
- 4 Choose Optimization from the list.
- 5 In the appearing window the optimization mode can be chosen from the list box as well as a delay time.

If you choose the Prefetch Sample Vial option, the next vial (for multiple injections the same vial) will be moved close to the injection port after the specified elapsed

Entering Settings

runtime. The injection cycle will not be started. This procedure reduces the total injection time without interfering with the current run.

By choosing the Overlap Injection Cycle option the next vial will be placed in the injection port after the selected elapsed runtime. The following restrictions apply when using the Overlap option:

- The elapsed runtime entry before the sample is processed must not interfere with the current running analysis.
- Overlapping works only for a given vial range (e.g. within one sequence line).
- The first run for a new vial range is done without overlap.
- Injector programs using valve switching commands must not be executed when overlapping is selected. They will fail with an error message. Since all injector programs created with the Agilent ChemStation require such commands in order to function correctly, none of them will work in the overlap mode.

The screen also allows you to enable/disable the option 'always keep transport arm next to last used vial'.

Setting up the Thermostatted Autosampler

If a thermostatted autosampler is part of your system you can enter specific setting using the Thermostat option in the settings screen.

Using this option you can set the temperature of the airflow into the autosampler. This setting is available in the Analysis screen, too. Activate the check box to have the temperature controlled as soon as the cooled autosampler is turned on.

Using the More ... option in the Settings screen and selecting Signals you can choose among several thermostat parameters (ambient temperature, heat sink, heat sink fan etc.) that will be referred to as "Autosampler: Auxiliary" in the Plot - Signals screen. This allows you to access a variety of thermostatted autosampler signals without overloading the Available Signals list box (see "Plot Screen" on page 30).

Aligning Transport Arm/Gripper

- 1 Press the Tests button [F3] in the System screen.
- 2 Choose Autosampler from the list.
- 3 Press the Align button [F1].
- 4 Choose Transport.
- 5 Enter the desired values in the corresponding fields.

The field X Correction sets the default value for the axis parallel to the front panel (left-right motion).

Theta Correction describes the default angle of the transport arm towards the front-back axis of the module (rotation).

For detailed information on the axes consult your Autosampler reference manual.

- 6 Select the Done button [F6] to accept settings.

The new values will be valid after performing the next hardware initialization or reset.

Tracking Autosampler History / EMF limits

- 1 Select the Records button [F4] in the System screen.
- 2 Select Autosampler from the list.
- 3 Press the Enter key.

You can check the number of 'Needle into Seat' injections the autosampler has made and the number of injection valve cycles. A limit can be specified for each function which you can use to schedule preventive maintenance.

Resetting the Autosampler

- 1 Select the Control button [F1] in the System screen.
- 2 Select Autosampler from the menu.
- 3 Press the Enter key.
- 4 Select the Reset button [F7] to reset the autosampler hardware.
This resets the injection valve, metering device and sampling unit.

Reset to Default Settings

- 1 Select the Settings button [F1] in the Analysis screen.
- 2 Select Autosampler from the menu.
- 3 Press the Enter key.
- 4 Select the Default button [F7] to reset the autosampler settings to their default values.

Troubleshooting the Autosampler

- 1 Select the Tests button [F3] in the System screen.
- 2 Select Autosampler from the menu.
- 3 Press the Enter key.

You can select a variety of injector steps to move individual parts of the autosampler.

The injector step functions can be used in troubleshooting to check the single steps of the injection cycle. For a complete description of each step, see the *Reference Manual* for the Agilent 1100 Series autosampler.

If you have a thermostatted autosampler, you can also test the auxiliary signals by pressing the Signal button in the Test screen. The auxiliary signals will be displayed graphically now (see “Setting up the Thermostatted Autosampler” on page 78).

Using the Manual Injection Valve

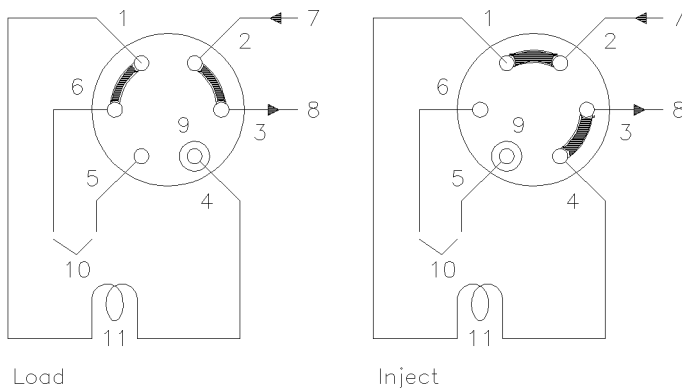
Operational details for the Agilent 1100 Series
manual injection valve

About the Injection Valve

The injection valve has six ports and can be moved manually between two positions. In the LOAD position, the mobile phase is led from the pump directly to the column, allowing you to inject your sample into the sample loop. In the INJECT position, the mobile phase is led through the sample loop, flushing the contents of the loop onto the column.

Figure 15

Valve Positions



Valve Positions:

- | | |
|---------------|--------------------|
| 1 to 6 | valve ports |
| 7 | inlet (from pump) |
| 8 | outlet (to column) |
| 9 | needle port |
| 10 | vents |
| 11 | sample loop |

Making An Injection

The sample loop can be either completely or partially filled. This depends on the amount of sample you have.

Completely Filling the Sample Loop

This is the conventional method in which an excess of sample is used to fill the sample loop completely. The volume of the loop determines the injection volume.

- 1 Fill syringe with sample.
 - 2 Move valve to LOAD position.
 - 3 Insert needle of syringe into needle port until needle touches stator face. Do not press too hard.
 - 4 Slowly inject sample.
 - 5 Leave syringe in position and move valve to INJECT position.
 - 6 Remove syringe.
-

Partially Filling the Sample Loop

Use this method when only small quantities of sample are available. In this method the syringe determines the injection volume.

- 1 With valve in INJECT position, use needle port cleaner to flush needle port with about 1 ml of mobile phase (this will reduce residual contamination from previous injection).
 - 2 Move valve to LOAD position.
 - 3 Fill syringe with required volume of sample (not more than half of loop volume).
 - 4 Insert needle of syringe into needle port until needle touches stator face. Do not press too hard.
-

How Much Sample Is Actually Injected?

- 5 Slowly inject sample.
- 6 Leave syringe in position and move valve to INJECT position.
- 7 Remove syringe.

How Much Sample Is Actually Injected?

When the sample loop is completely filled, the amount of sample injected is equal to the volume of the sample loop plus the volume of the valve passages (1 in rotor and 2 in stator). This means that the actual amount of sample injected will be different to the nominal value designated to your sample loop. However, since both standards and samples are analyzed using the same loop, you rarely need to know the absolute volume of the loop.

If you do need to know the actual volume of a sample loop, we recommend you calibrate it fitted to the valve, so that you also take the valve passages into account.

How Much Sample Do I Need?

To completely fill the sample loop an excess of sample is required. This is about 2 to 3 loop volumes of sample to achieve 95 % of the maximum loop volume (the remainder is residual mobile phase in the sample loop). Determine the optimum number of loop volumes experimentally for your particular application.

When you partially fill the sample loop, do not inject more than half of the sample volume. As you inject your sample into the loop, the sample mixes with the solvent already in the loop and some of the sample could be lost through port 6 if you try to inject too much.

An Alternative Way to Fill The Loop

When the sample loop is completely filled you can either inject the sample into the loop or use the syringe to draw the sample through the loop.

- 1** Move valve to LOAD position.
- 2** Place vent tube from port 6 into sample vial.
- 3** Insert needle of syringe into needle port until needle touches stator face. Do not press too hard.
- 4** Slowly draw sample into syringe.
- 5** Leave syringe in position and move valve to INJECT position.
- 6** Remove syringe.

If you use this method, always flush the loading passages (vent tube on port 6) after each injection to prevent cross-contamination between injections. Remember — to flush port 6 and the vent tube, the valve must be in the LOAD position.

Using the Manual Injection Valve
An Alternative Way to Fill The Loop

Using the Detectors

Operational details for the Agilent 1100 Series
variable wavelength, multiple wavelength, refractive
index, fluorescence light, and diode array detectors

Turning the Lamp On and Off

To turn the lamp on and off use the On/Off button [F7] in the Analysis screen:

- 1 Enter the Analysis screen.
- 2 Select the On/Off button [F7].
- 3 Select the Lamp button [F6].

The same can be achieved by using the On/Off button [F8] in the Detector Settings or Control screen available through the Analysis or System screens.

NOTE

When the lamp is turned on, a lamp ignition routine is started. For more information see “Lamp Ignition Routine (VWD only)” on page 92.

Entering Settings

All detector settings can be entered in the Detector Settings screen. You can access this screen by selecting the Settings button [F1] in the Analysis screen and choosing the appropriate detector. Commonly used settings such as wavelength can be set directly in the Analysis screen.

Diode Array Detector

The wavelength for signal A can be entered in the Analysis screen. The sample wavelength, reference wavelength, bandwidth, peakwidth and slit width can be entered in the DA Detector Settings screen. This screen is accessed by selecting the Settings button [F1] in the Analysis screen. Additional entries (autobalance etc.) and the spectrum settings (range, threshold) are accessed by selecting the More... button [F1] under DA Detector Settings.

Variable Wavelength Detector

Settings for this detector include the wavelength, peakwidth and signal polarity. Pressing the More ... button [F1] and choosing Settings you can enter autobalance

settings and the margin for negative absorbance. The option Spectrum allows you to enter a wavelength range for a sample spectrum.

Fluorescence Light Detector

For the fluorescence light detector the following settings can be entered: Excitation wavelength, emission wavelength and Multi wavelength settings can be set in the regular Settings screen. With the More ... button [F1] you can specify further detector settings, i.e. peakwidth and PMT-Gain (baseline behaviour, reference, polarity, spectral range fit), multi wavelength settings (spectrum, excitation range), 3D scanning settings (excitation and emission ranges) and phosphorescence detection mode settings.

Multiple Wavelength Detector

The wavelength for signal A can be entered in the Analysis screen. The wavelength for signal B, slit width and peak width can be entered in the MW Detector Settings screen. This screen is accessed by selecting the Settings button [F1] in the Analysis screen. Additional signals and functions are accessed by selecting the More... button [F1] under MW Settings.

NOTE

Only the G1323B version of the control module allows control of the Multiple Wavelength Detector. The G1323A will display "resident or unsupported module" if a RI detector is configured in the Agilent 1100 system.

Refractive Index Detector

The control module allows to set the RID cell temperature in the Analysis screen.

Temperature, Peakwidth, polarity and automatic recycling can be set in the RI detector settings screen. This screen is accessed by selecting the Settings button [F1] in the Analysis screen.

Automatic zero and automatic purge can be set by selecting the More... button [F1] under RI Settings.

NOTE

Only the G1323B version of the control module allows control of the Refractive Index Detector. The G1323A will display "resident or unsupported module" if a RI detector is configured in the Agilent 1100 system.

Resetting the Baseline

You can reset the baseline using the Balance button [F1] (Zero button [F1] when using the Refractive Index Detector) in the Detector Control screen. Please note that this procedure is not available for the Fluorescence Light Detector:

- 1 Select the Control button [F1] in the System screen.
- 2 Select the Detector from the menu.
- 3 Press the Enter key.
- 4 Select the Balance button [F1].

You can achieve the same using the m (menu) key in the Detector Settings window available from the Analysis screen.

Configuring the Detector

- 1 Select the Configure button [F2] in the System screen.
- 2 Select the Detector from the list.
- 3 Press the Enter key.

You can now configure the detector analog output(s), lamp-on at power-on and the interfaces.

Troubleshooting the Detector

- 1 Select the Tests button [F3] in the System screen.
- 2 Select the Detector from the menu.
- 3 Press the Enter key.
- 4 Select the test you require.

The selection of tests depends on the. For information about each test, see the *Reference Manual* for the Agilent 1100 detectors.

NOTE

The full test capability is only available from the LC ChemStation.

Tracking Detector History

- 1 Select the Records button [F4] in the System screen.
- 2 Select the Detector from the menu.
- 3 Press the Enter key to display the logbook.

You can check, change and reset the lamp burn time and number of ignitions. For the FL detector the flash lamp lifetime and for the RI detector the time since last purge can be checked.

Resetting the Detector

- 1 Select the Settings button [F1] in the Analysis screen.
 - 2 Select the Detector from the menu.
 - 3 Press the Enter key.
 - 4 Select the Default button [F7] to reset the detector settings to their default values.
-

Lamp Ignition Routine (VWD only)

When the lamp is turned on, the following routine is implemented:

- 1-minute warm-up, within this time the grating position is re-initialized.
Check of wavelength setting at 656 nm emission (must be within 1 nm) at reference side only (to have no influence from flow cell condition).
If not correct, a message for calibration is displayed. Perform a wavelength calibration (available within VWD Tests).
 - Check of intensity
 - at 250 nm without cut-off filter inserted,
 - at 250 nm with cut-off filter inserted.If the intensity check is not OK, the cut-off filter operation is not OK.
-

Configuring the Analog Channel Output

You may use the LC Series 1100 detectors to test or monitor optical parameters on external analog equipment connected via the analog output (Fluorescence Light, Diode Array and Multiple Wavelength Detectors have two analog outputs, Variable

Configuring the Analog Channel Output

Wavelength Detector and Refractive Index Detector have one) at the back side of the module.

To configure the analog outputs:

- 1** Press the Settings button [F1] in the Analysis screen.
- 2** Choose the detector and press enter.
- 3** Press the Analog button [F3].
- 4** Either select the signal you want to monitor from the signal source drop down list. Or, if you have a variable wavelength detector (fluorescence light detector), choose the signal(s) from the list(s) to be routed to the analog output(s).
- 5** Enter the values for Zero Offset (sets a baseline offset in order to recognize negative drifts) and Attenuation (sets an absorbance range) for each of the two connectors.
- 6** Press the Done button [F6].

Using the Detectors

Configuring the Analog Channel Output

Using the Column Compartment

Operational details for the Agilent 1100 Series
thermostatted column compartment

Turning the Column Compartment On and Off

You can turn the column compartment on and off by using the On/Off button [F7] in the Analysis screen:

- 1 Enter the Analysis screen.
- 2 Select the On/Off button [F7].
- 3 Select the Temp button [F7].

The same can be achieved through the Settings or Control buttons [F1] in the Analysis or Systems screens.

Entering Settings

The temperature can be entered in the Analysis screen. The temperature (for left and/or right heater) can be entered in the Column Compartment Settings screen. This screen is accessed by selecting the Settings button [F1] in the Analysis screen. Additional settings are accessed by selecting the More... button [F1] under Column Compartment Settings.

Configuring the Column Compartment

- 1 Select the Configure button [F2] in the System screen.
- 2 Select Column Comp from the list.
- 3 Press the Enter key.

You can now configure the leak-detection mode, temperature-on at power-on, the interfaces and the column identification module.

Configuring the Column ID Module

- 1 Select the Configure button [F2] in the System screen.
- 2 Select Column Comp from the list.
- 3 Press the Enter key to enter the Column Compartment Configuration screen.
- 4 Press the Column ID button [F4] to enter Column ID screen.
- 5 To toggle between left and right column press the Left/Right button [F8].

NOTE

If no column tag is sensed, the record fields are not active and the left/right tag sign in the upper right corner of the window is crossed out.

- 6 Enter your column data into the fields as required. Additional fields are available by pressing the More ... button [F1].
- 7 Pressing the Write button [F7] will transfer the information into the column tag.
- 8 Pressing the Right button [F8] will show the information fields of the right column tag (if column is installed).

NOTE

The column ID information is updated as soon as a column with ID tag is installed correctly. So you can check the number of injections that have been made on the column in this screen.

Selecting Separated or Combined Mode

This mode allows the temperature setting of both heaters independent from each other. If not enabled both heaters are kept on the same temperature.

- 1 Select the Settings button [F1] in the Analysis screen.
- 2 Select Column Comp from the menu.
- 3 Enable the Separated mode by selecting the check box.
- 4 To enable the Combined mode, deselect the check box.

Selecting the Column Switching Valve (optional)

The column switching valve is optional.

- 1 Select the Settings button [F1] in the Analysis screen.
- 2 Select Column Comp from the menu.
- 3 Press the Column Switch button.
- 4 Select mode 1 or 2.
- 5 Press the Done button [F6] and leave this screen.

Troubleshooting the Column Compartment

- 1 Select the Tests button [F3] in the System screen.
- 2 Select Column Comp from the menu.
- 3 Press the Enter key.
- 4 Select the test you require.

You can select the Calibrate test to check the operation of the Column Compartment. For information about the test, see the *Reference Manual* for the Agilent 1100 Series column compartment.

Tracking Column Compartment History

- 1 Select the Records button [F2] in the System screen.
 - 2 Select Col. Comp from the list.
 - 3 Press the System or Maint. Log button [F4] to display the logbook.
-

Tracking Column ID History

- 1 Select the Configure button [F4] in the System screen.
- 2 Select Column Comp from the list.
- 3 Press the Enter key to enter the Column Compartment Configuration screen.
- 4 Press the Column ID button [F2] to enter the Column ID screen.

NOTE

If no column tag is sensed, the parameter entry fields are not active, the left/right tag sign is crossed out and the More ... button is unavailable.

Resetting the Column Compartment

By selecting the More ... button [F1] you can enter values and check a variety of features for the installed column. For example, the maximum pressure allowed and the maximum recommended temperature.

You can check the number of injections that have been made on the column.

Resetting the Column Compartment

- 1 Select the Settings button [F1] in the Analysis screen.
- 2 Select Column Comp from the menu.
- 3 Press the Enter key.
- 4 Select the Default button [F7] to reset the detector settings to their default values.

Part 3

9	Running an Isocratic Analysis	103
10	Running a Gradient Analysis	109
11	Running Multiple-Vial Analyses	117
12	Running an Injector Program	129

Using the Agilent 1100 Series LC System With Control Module

Running an Isocratic Analysis

How to analyze the Agilent Technologies isocratic standard sample using a single injection analysis

What You Will Need

Instruments	Agilent 1100 Series isocratic, binary or quaternary pump, autosampler and a UV-detector.
Column	A 125 mm × 4.0 mm Hypersil ODS, 5 µm (Agilent Technologies part number 7982618-564).
Solvents	For the isocratic pump, a solvent mixture of LC grade bidistilled water (35 %) and acetonitrile (65 %).
Sample	The Agilent Technologies isocratic standard sample (Agilent Technologies part number 01080-68704). This contains 0.15 wt.% dimethylphthalate, 0.15 wt.% diethylphthalate, 0.01 wt.% biphenyl and 0.03 wt.% o-terphenyl dissolved in methanol.

Preparing the LC System

- 1 For the isocratic pump, fill the solvent bottle with the mixture of LC-grade bidistilled water (35 %) and acetonitrile (65 %). For the binary or quaternary pump, fill one solvent bottle with bidistilled water (channel A) and the other with acetonitrile (channel B).
- 2 Turn on the detector lamp and pump using the On/Off button [F7] in the Analysis screen. (Use Action keys then to select the module)
- 3 For the quaternary pump, turn on the degasser by pressing the line-power switch.
- 4 Purge the pump. For more information see Chapter 3 “Using the Pump”.
- 5 Allow the detector at least 15 minutes to provide a stable baseline.
- 6 Fill the contents of a Agilent Technologies isocratic standard sample ampoule into a vial and seal the vial with a cap. Place the vial in position 1 of the autosampler tray.
- 7 Pump the water/acetonitrile (35/65 %) mobile phase through the column for 10 minutes at a flow rate of 2 ml/min.

Entering Settings

To set up the isocratic analysis you will set the LC system settings to default and then modify selected settings, the other settings will remain with their default values. You will then save these settings to a method called ISO.

- 1 Enter the Analysis screen.
- 2 Set the vial range as 1 to 1.
- 3 Set the number of injections to 1.
- 4 Press Enter to skip the method name section.
- 5 Select the Settings button [F1].
- 6 Select LC System from the menu.
- 7 Select the Default button [F7] and select the Yes button to load defaults.
- 8 Press the Esc key.
- 9 Select the Settings button [F1].
- 10 Select the pump from the menu (either ISO, BIN or QUAT).
- 11 Press the Pressure button [F3].
- 12 Enter 400 as the upper pressure limit.
- 13 Select the Done button [F6].
- 14 Enter the following values: %B 65, (%C OFF, %D OFF for quaternary pump. If you have a binary pump, set %B to 65.).
- 15 Set the Flow to 1,5 ml/min.
- 16 Confirm by selecting “Done” to enter the Analysis screen
- 17 Enter the values shown in Table 2 in the Analysis screen.

NOTE

The channels of a pumping system are named A, B, C and D (depending on the pump type). %A is automatically calculated by $100\% - (\%B + \%C + \%D)$. If no values for %B, %C and %D are entered, %A is always 100%.

Table 2

Values in Analysis Screen	
Setting	Value
Detection wavelength	254 nm
Injection volume	1 µl
Stoptime	6 min

Saving Settings in a Method

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select the Save As button [F8].
- 3 Enter the method name as ISO using the selection keys (also see “Specifying a Method Name” on page 39)
- 4 Press the Done button [F6] to save the method.
- 5 Press the Esc key to return to the Analysis screen.

Observing the Chromatogram

- 1 Select the Plot button [F6] in the Analysis screen.
- 2 Press the Select button [F6].
- 3 Choose a Signal from the Available Signals list box.
- 4 Press the Enter key

NOTE

You can choose several signals at a time. The plot function will display all signals that are shown in the Selected Signals list box. Use the selection keys to navigate within the list boxes and press the Enter key to move Signals from one box to the other.

- 5 Select the Done button [F6] to display the chromatogram.
- 6 Press the Esc key to go back to the Analysis screen.
- 7 Press the Start button [F8].
- 8 Select OK to confirm the vial range and injection number and press the Enter key to start the analysis.
- 9 Press the Plot button [F6] to show the chromatogram

A typical chromatogram for this analysis is shown in Figure 16.

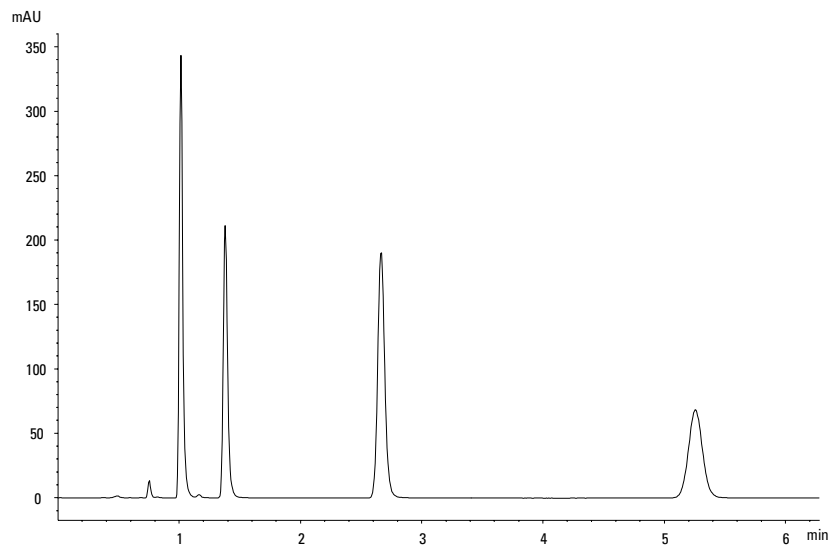
The exact profile of the chromatogram will depend on the column you have used. Differences in retention times and areas of the peaks in your chromatogram and the one shown in Figure 16 might be a result of variations in the concentration of the sample from batch to batch, the quality of the solvents used and the column temperature.

NOTE

You can rescale the plot using the Rescale button [F7], or the cursor keys or you define the plot window within the Setup.

Figure 16

Analysis of Isocratic Standard Sample



Starting the Analysis

NOTE

The analysis can be started from various other screens, e.g. System, Status, Sampels or Sequence screen.

- 1 Enter the Analysis screen.
- 2 Select the Start button [F8].
- 3 Select OK to confirm the vial range and injection number and press the Enter key to start the analysis.

Running a Gradient Analysis

How to analyze the Agilent Technologies isocratic standard sample using a gradient analysis

What You Will Need

Instruments	Agilent 1100 Series binary or quaternary pump, autosampler, UV-detector and degasser.
Column	A 125 mm × 4.0 mm Hypersil ODS, 5 μm (Agilent Technologies part number 7982618-564).
Solvents	LC-grade bidistilled water and acetonitrile.
Sample	The Agilent Technologies isocratic standard sample (Agilent Technologies part number 01080-68704). This contains 0.15 wt.% dimethylphthalate, 0.15 wt.% diethylphthalate, 0.01 wt.% biphenyl and 0.03 wt.% o-terphenyl dissolved in methanol.

Preparing the LC System

- 1 Fill one solvent bottle with bidistilled water (channel A) and the other with acetonitrile (channel B).
- 2 Turn on the detector lamp and pump using the On/Off button [F7] in the Analysis screen.
- 3 For the quaternary pump, turn on the degasser by pressing the line-power switch.
- 4 Purge the pump. For more information see Chapter 3 “Using the Pump”.
- 5 Allow the detector at least 15 minutes to provide a stable baseline.
- 6 Fill the contents of a Agilent Technologies isocratic standard sample ampoule into a vial and seal the vial with a cap. Place the vial in position 1 of the autosampler tray.
- 7 Pump the water/acetonitrile (35/65 %) mobile phase through the column for 10 minutes at a flow rate of 2 ml/min.

Entering Settings

To set up the gradient analysis you will set the LC system settings to default and then modify selected settings, the other settings will remain with their default values. You will then save these settings to a method called GRAD.

To set up the solvent gradient, you will create a timetable in the method. This is done using the Timetable screen, accessed by selecting the Timetable button [F2] in the Pump settings screen.

The example assumes that the Timetable is empty. If the Timetable is not empty, use the All Lines item, accessed by pressing the Delete button.

Entering Settings

- 1 Enter the Analysis screen.
- 2 Set the Vial Range as 1 to 1.
- 3 Set the Number of Injections to 1.
- 4 Skip the method name field
- 5 Select the Settings button [F1].
- 6 Select LC System from the menu.
- 7 Press the Enter key.
- 8 Select the Default button [F7] and select the Yes button to load defaults.
- 9 Press the Esc key to enter the Analysis screen.
- 10 Enter the following values in the Analysis screen.

Table 3**Values in the Analysis Screen**

Setting	Value
Flow	1.5 ml/minute
Detection wavelength	254 nm
Injection volume	1 µl
Stoptime	5 min

- 11 Select the Settings button [F1].
- 12 Select BIN or QUAT pump from the menu.
If you have a quaternary pump set %B to 65, %C and %D to OFF.
If you have a binary pump set %B to 65.

NOTE

The channels of a pumping system are named A, B, C and D (depending on the pump type). %A is automatically calculated by $100\% - (\%B + \%C + \%D)$. If no values for %B, %C and %D are entered, %A is always 100%.

Entering Settings

- 13 Select the Pressure button.
- 14 Set the upper pressure limit to 400 bar.
- 15 Select the Done button.
- 16 Select the Timetable button.
- 17 Select the Insert button.
- 18 Enter a value of 2 minutes and select Composition as the setting.
If you have a quaternary pump, set the %B composition to 65, %C and %D to 0.00.
If you have a binary pump, set the %B composition to 65.
- 19 Select the Enter button to accept entries (Action Key F7).
- 20 Enter the following information for another two lines.
time 4 : %B = 95
time 5 : %B = 65
- 21 Press the Done button and verify your timetable entries.
- 22 Press the Esc key until you are in the Analysis screen.

NOTE

Use Enter to store timetable lines.
Use Done when all lines have been entered.
By pressing the Esc key you return to the previous screen.

Saving Settings in a Method

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select the Save As button [F8].
- 3 Enter the method name as GRAD using the selection keys (also see “Specifying a Method Name” on page 39)
- 4 Press the Done button [F6] to save the method.
- 5 Press ESC to return to the Analysis screen.

Observing the Chromatogram

- 1 Select the Plot button [F6] in the Analysis screen.
- 2 Press the Select button [F6].
- 3 Choose a Signal from the Available Signals list box.
- 4 Press the Enter key

NOTE

You can choose several signals at a time. The plot function will display all signals that are shown in the Selected Signals list box. Use the selection keys to navigate within the list boxes and press the Enter key to move Signals from one box to the other.

- 5 Select the Done button [F6] to display the chromatogram.
- 6 Press Esc to go back to the Analysis screen.
- 7 Press the Start button [F8].
- 8 Select OK to confirm the vial range and injection number and press the Enter key to start the analysis.
- 9 Press the Plot button [F6] to show the chromatogram.

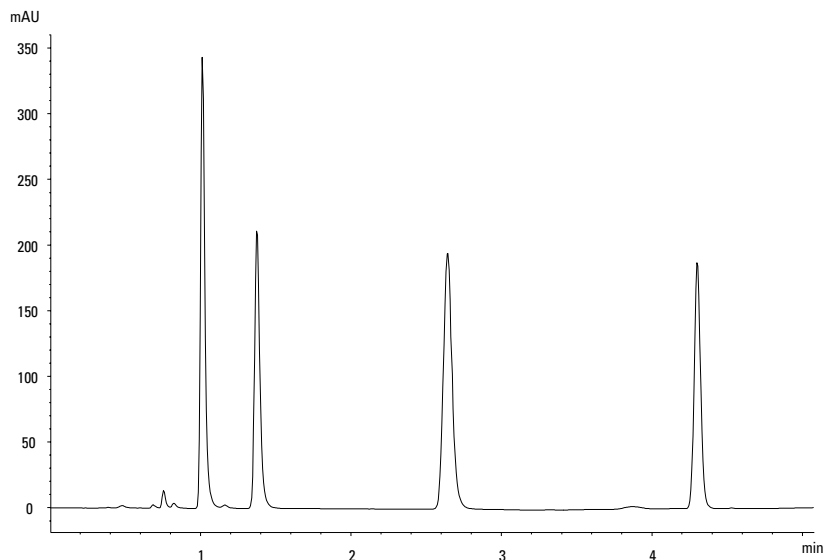
Running a Gradient Analysis

Observing the Chromatogram

A typical gradient chromatogram for this analysis is shown in Figure 17.

Figure 17

Gradient Analysis of Isocratic Standard



The exact profile of the chromatogram will depend on the column you have used. Differences in retention times and areas of the peaks in your chromatogram and the one shown in Figure 17 might be a result of variations in the concentration of the sample from batch to batch, the quality of the solvents used and the column temperature.

If you compare this chromatogram with the one from Chapter 9 “Running an Isocratic Analysis” you will notice the solvent gradient has reduced the elution time of the fourth peak in the chromatogram.

NOTE

You can rescale the plot using the Rescale button [F7], or the cursor keys or you define the plot window within the Setup.

Starting the Analysis

NOTE

The analysis can be started by pressing the Start button [F8] from the Samples screen, the System screen, or the Status screen.

- 1 Enter the Analysis screen.
- 2 Select the Start button [F8].
- 3 Select OK to confirm the vial range and injection number and press the Enter key to start the analysis.

Running Multiple-Vial Analyses

How to setup multiple vial analyses using the same method and different methods

Analyzing Multiple Vials Using the Same Method

This section describes how to set up a 25-vial analysis with one injection from each vial. You will use a previously created method. The samples are located in positions 1 to 25 of the autosampler tray. For details see “Working with Methods” on page 38.

- 1 Enter the Analysis screen.
- 2 Set the Vial Range from 1 to 25.
- 3 Set the number of injections to 1.
- 4 Select the method you want to use with the scroll buttons (▲ ▼).
- 5 Select the Start button [F8].
- 6 Select OK to confirm the vial range and injection number and press the Enter key to start the analysis.

Analyzing Multiple Vials Using Different Methods

This section describes how to set up a 50-vial analysis using three methods which you have previously created called e.g. METH1, METH2 and METH3. For example: METH1 and METH2 have the same analytical settings but differ in the injection volume and stoptime values. METH3 uses a different temperature and requires a wait time of 30 minutes for the LC system to equilibrate.

NOTE

This way of combining methods is called *Sequence*.

- The first 20 vials are analyzed using METH1 with one injection per vial,
- the next 20 vials are analyzed using METH2 with two injections per vial,
- the last 10 vials are analyzed using METH3 with three injections per vial.

The vials are located in positions 1 to 50 of the autosampler tray.

Running Multiple-Vial Analyses

Analyzing Multiple Vials Using Different Methods

The example assumes that the sequence table is empty. If the sequence table is not empty, use the All Lines item, accessed by pressing the Delete button [F6].

- 1 Select the Sequence [F4] in the Analysis screen.
- 2 For sequence line 1, select the Insert button [F7] and enter:

Vial Range	1 to 20
#Inj. 1	1
Inj. Vol.	Default
Method	Meth1

- 3 Select the Enter button to accept entries.

- 4 For sequence line 2, enter:

Vial Range	21 to 40
#Inj.	2
Inj. Volume	Default
Method	METH2

- 5 Select the Enter button to accept entries.

- 6 For sequence line 3, enter:

Vial Range	41 to 50
#Inj.	3
Inj. Volume	Default
Method	METH3
Wait	30 minutes

- 7 Select the Done button [F6] to enter sequence line 3 settings and complete the sequence.
- 8 Select the Start button [F8] in the Sequence screen.
- 9 Select the OK button to start the sequence.

Single-Level Calibration Sequences

The following procedure describes how to set up a calibration sequence for an analysis which uses single-level calibration.

There is one calibration standard (C) and 9 samples (S).

The analysis requires that:

- each sample is analyzed in duplicate,
- the calibration standard is analyzed once before the samples and re-analyzed once after every 2 samples,
C S S C S S C S S C S S C S C
- the calibration standard is located in position 90 and the 9 sample vials are in positions 1 to 9 of the autosampler tray, and
- the method called METH1 is used for the samples and standards.

The example assumes that the sequence table is empty. If the sequence table is not empty, use the All Lines item, accessed by pressing the Delete button [F6].

- 1 Select the Sequence button [F4] in the Analysis screen.
- 2 For sequence line 1, select the Insert button [F7] and enter:

Vial Range	1 to 9
#Inj.	2
Inj. Volume	Default
Method	METH1

Running Multiple-Vial Analyses

Multiple-Level Calibration Sequences

- 3 Select the Calibration button [F1] to display the Calibration Settings screen for sequence line 1 and choose Edit.
- 4 Enter the following information:

Calibration vial range	90 to 90
Number of injections	1
Calibration method	METH1
Recalibrate every	2 vials multi
Before	On
After	On
- 5 Press the Done button [F6] until the Sequence screen appears.
- 6 Select the Start button [F8] in the Sequence screen.
- 7 Press the Enter key to start the sequence.

Multiple-Level Calibration Sequences

The following sections describe how to set up calibration sequences for analyses which use multiple-level calibration.

Recalibrating With the Same Group of Standards

There are three calibration standards of different concentrations (C1, C2, C3) and 15 samples (S). The standards and samples are analyzed using the same method.

The analysis requires that:

Running Multiple-Vial Analyses

Recalibrating With the Same Group of Standards

- each sample is analyzed once,
- the calibration standards are analyzed twice before the samples and re-analyzed twice after every 5 samples,
C1 C2 C3 S10-S14 C1 C2 C3 S15-S19 C1 C2 C3 S20-S24 C1 C2 C3
- the calibration standards are located in positions 90 to 92 of the autosampler tray,
- the 15 sample vials are located in positions 10 to 24 of the autosampler tray, and
- the samples and standard are analyzed using a method called METH1.

The example assumes that the sequence table is empty. If the sequence table is not empty, use the All Lines item, accessed by pressing the Delete button [F6].

- 1 Select the Sequence button [F4] in the Analysis screen.
- 2 For sequence line 1, select the Insert button [F7] and enter:

Vial Range	10 to 24
#Inj	2
Inj. Volume	Default
Method	METH1
- 3 Select the Calibration button [F1] to display the Calibration Settings screen for sequence line 1 and choose Edit.
- 4 Enter the following information:

Calibration vial range	90 to 92
Number of injections	2
Calibration method	METH1
Recalibrate every	5 vials multi
Before	On
After	Off
- 5 Select the Done button [F6] to accept entries.
- 6 Select the Start button [F8] in the Sequence screen.
- 7 Press the Enter key start the sequence.

The autosampler now analyzes:

- the three calibration standards in duplicate,
- sample vials 10 through 14,
- the three calibration standards in duplicate,
- sample vials 15 through 19,
- the three calibration standards in duplicate,
- sample vials 20 through 24, and
- the three calibration standards in duplicate.

Recalibrating With Multiple Groups of Standards

There are two different types of samples, A and B that need to be analyzed.

The analysis for sample type A requires a 5 µl injection and a stoptime of 8 minutes.

The analysis of sample type B requires a 2 µl injection and a stoptime of 5 minutes.

For sample type A:

- there are 3 calibration standards of different concentrations and 6 samples,
- each sample must be analyzed once,
- the calibration standards must be analyzed in duplicate and re-analyzed after every 2 samples,
C1 C2 C3 S7 S8 C1 C2 C3 S9 S10 C1 C2 C3 S11 S12 C1 C2 C3
- The calibration standards of type A are in positions 1, 2 and 3 of the autosampler tray and the 6 sample vials are in positions 7 to 12, and
- the samples and the calibration standards use the same method called METH1.

For sample type B:

- there are 3 calibration standards of different concentrations and 9 samples,
- each sample must be analyzed once,
- the calibration standards must be analyzed twice and re-analyzed after every 3 samples,

Recalibrating With Multiple Groups of Standards

C1 C2 C3 S13-S15 C1 C2 C3 S16-S18 C1 C2 C3 S19-S21 C1 C2 C3

- the calibration standards of type B are in positions 4, 5 and 6 of the autosampler tray and the 9 sample vials are in positions 13 to 21, and
- the samples and calibration standards of type B use different methods.

The samples use METH2 and the calibration standards use METH3. These methods contain the same analytical parameters and differ only in the analysis stoptime.

The example assumes that the sequence table is empty. If the sequence table is not empty, use the All Lines item, accessed by pressing the Delete button [F6].

1 Select the Sequence button [F4] in the Analysis screen.

2 For sequence line 1, select the Insert button [F7] and enter:

Vial Range	7 to 12
#Inj.	1
Inj. Volume	Default
Method	METH1

3 Select the Calibration button [F1] to display the Calibration Settings screen for sequence line 1 and choose Edit.

4 Enter the following information:

Calibr. Vial range	1 to 3
#Inj.	2
Inj. Volume	Default
Calibr. method	METH1
Recalibrate every	2 vials multi
Before	On
After	Off

5 Select the Done button [F6] to accept entries.

6 Move the highlighted bar to line 2 (by pressing the Selection key down) and press the Enter key.

7 For sequence line 2 enter:

Vial Range	13 to 21
#Inj.	1

Recalibrating With Multiple Groups of Standards

Inj. Volume Default

Method METH2

- 8 Select the Calibration button [F1] to display the Calibration Settings screen for sequence line 2 and choose Edit.

- 9 Enter the following information:

Calibr. vial range 4 to 6

#Inj 2

Inj. Volume Default

Calibr. method METH3

Recalibrate every 3 vials multi

Before On

After Off

- 10 Select the Done button [F6] to accept entries.

- 11 Select the Start button [F8] in the Sequence screen.

- 12 Press the Enter key to start the sequence.

The autosampler now analyzes:

- three type A calibration standards in duplicate,
- type A samples in vials 7 and 8,
- three type A calibration standards in duplicate,
- type A samples in vials 9 and 10,
- three type A calibration standards in duplicate,
- type A samples in vials 11 to 12,
- three type A calibration standards in duplicate,
- three type B calibration standards in duplicate,
- type B samples in vials 13, 14 and 15,
- three type B calibration standards in duplicate,
- type B samples in vials 16,17 and 18,
- three type B calibration standards in duplicate,
- type B samples in vials 19, 20 and 21, and

- three type B calibration standards in duplicate.

Synchronizing Analyses with External Devices

With an APG remote connector the system can be connected to external devices in order to synchronize the analyses. This is necessary when an external device needs some time in order to get ready for a new analysis and when transmitting of a start request is required (see “Connecting External Devices” on page 51 for further information on interfaces).

When executing an analysis by pressing the Start button [F8] you will see a window where sequence modes can be chosen.

In any case, do all the analysis preparation using the control module.

NOTE

A “Start” command is used to start the analytical run from the point of injection and is usually issued by the autosampler.

A “Start Request” command causes the autosampler to take the next vial and place it under the injection needle (also see “APG Remote” on page 51 and “Optimizing Autosampler Performance” on page 77)

The “Start button” [F8] on the control module is used to start a vial range or sequence analysis.

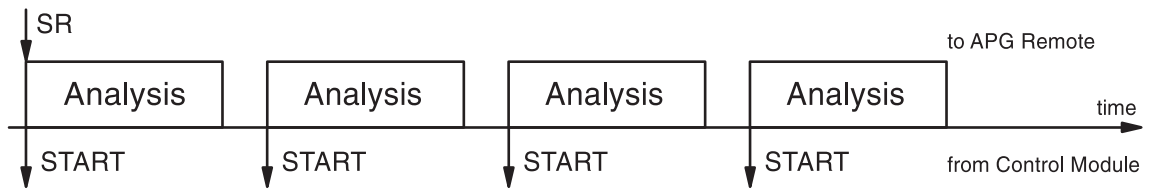
Standard

In the standard mode the analysis is under the command of the control module. The control module will issue a Start Request command to the autosampler as soon as all modules are ready for the next analysis. The autosampler issues the Start command at the point of injection. With an Agilent 1100 Autosampler integrated in the system and no external devices this is the normal operation mode.

Send Single Start Request

After you start the analysis with the control module it will generate a single start request on the APG remote lines. This allows to trigger the external device which will start each injection by sending a start signal. The vial range or sequence will be started by the control module, but for every new injection the external device must give the start command.

Figure 18 **Send single external start request**

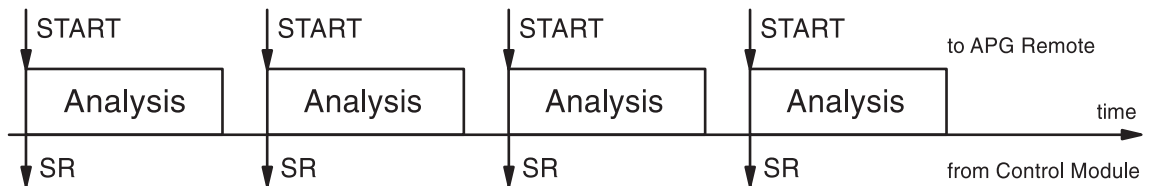


In this mode the control module will simply track the progress of the vial range or sequence. It will indicate the proper sample, change sequence lines, load methods or other sequence related actions for each run.

Send Repeated Start Request

This will cause the control module to generate start requests before each run. The external device starts each injection then by sending a start signal to the APG remote line. That is, after all the programming has been completed on the control module and the Start button [F8] has been pressed, a Start Request is issued before each run and the external device must give the Start command for the injection process.

Figure 19 **Send repeated external start request**

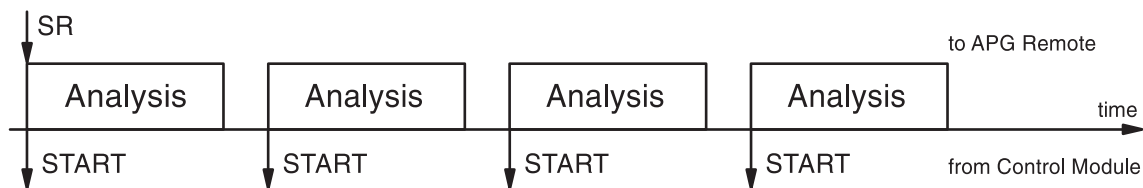


Here, too, the module will simply track the progress of the analysis.

Wait for Single (External) Start Request

After pressing the Start button [F8] the autosampler will wait for a single external start request on the APG remote lines. When the start request is received, the complete vial range or sequence is done as in standard mode under the command of the control module.

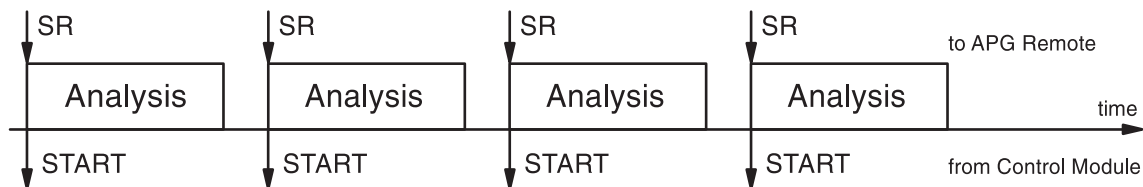
Figure 20 Wait for single external start request



Wait for Repeated Start Request

After pressing the Start button [F8] the autosampler will wait for external start requests before each vial in the range or sequence. They have to be generated by the external device. This mode is necessary when the external device needs extra time to get ready for the next analysis and thus has to be in charge of the start event.

Figure 21 Wait for repeated external start request



NOTE

An Agilent 1100 variable wavelength detector or diode array detector will perform a balance (provided that Auto Balance is set to Prerun in the More ... Settings screen) when receiving a start command from the control module. This will only happen in the Standard and Wait for single (repeated) start request modes. In the Send single (repeated) start request modes a balance before the run will NOT be performed. If regular balancing is required, set the Auto Balance check box to Postrun.

Running an Injector Program

How to create an injector program

Creating an Injector Program

The injector program is part of the method. The injector program screen can be accessed using the Inj. Program button [F3] in the Autosampler Settings screen.

This section describes an injector program to do sample preparation involving a precolumn derivatization reaction. This is required when the analytes lack chromophores and in their original chemical structure cannot be detected with the required sensitivity.

To create this injector program you will:

- set the LC system to the default settings
- modify the injector program settings in the Inj. Program screen, the other settings will remain with their default values, and
- save the method with the name DERIV.

When the method called DERIV is started, the program is executed followed by the analytical run.

The first step of the program involves drawing derivatization reagent into the needle capillary, followed by sample and then reagent again. To ensure the reagent and sample react efficiently, the plunger in the metering device of the autosampler is moved back and forth. This mixes the sample and reagent as they travel up and down the capillary. Following mixing, the derivatization requires a pause to allow the sample and reagent to fully react, before the derivatized sample is injected onto the column.

Entering Injector Program Settings

- 1 Enter the Analysis screen.
- 2 Select the Settings button [F1].
- 3 Select LC System from the menu.
- 4 Select the Default button [F6] and select the Yes button to load defaults.
- 5 Press Esc to go back to the Analysis menu.

Saving the Method

- 6 Select the Settings button [F1].
- 7 Select Autosampler from the menu.
- 8 Select the Inj.Program button [F3] to display the Program screen.
- 9 For program line 1, select the Insert button [F7].
- 10 Select the DRAW function and enter the draw settings as described in Table 4.
- 11 Select the Enter button [F7] to accept entries.
- 12 Repeat the procedure for the other functions to complete the program.
- 13 Press the Esc key until you are back in the Settings screen.

Injector Program Lines

Table 4 on page 132 describes the program lines which will complete the required derivatization reaction.

Saving the Method

- 1 Select the Method button [F3] in the Analysis screen.
- 2 Select the Save As button [F8].
- 3 Enter the method name as DERIV using the selection keys (also see “Specifying a Method Name” on page 39)
- 4 Press the Done button [F6] to save the method.
- 5 Press the Esc key to return to the Analysis screen.

The injector program is now part of the method called DERIV and can be started by selecting the Start button [F8] in the Analysis screen.

Table 4

Injector Program Lines

Line	Function	Parameters	Comment
1	DRAW	2 μ l (AMOUNT) from Air (SOURCE)	Draw 2 μ l from air into the needle capillary, to separate the mobile phase sitting in the capillary from the sample.
2	DRAW	1 μ l (AMOUNT) from vial 1(SOURCE)	Draw 1 μ l from derivatization reagent from vial #1 in the autosampler tray into the needle capillary.
3	WASH	in vial 2 (VIAL), 1 time (Cycles)	Wash the needle tip in the wash vial #2.
4	DRAW	2 μ l (AMOUNT) from sample (SOURCE)	Add 2 μ l of the sample in the autosampler tray to the derivatization volume already in the needle capillary.
5	WASH	in vial 2 (VIAL), 1 time (CYCLE)	Wash the needle tip in the wash vial #2.
6	DRAW	1 μ l (AMOUNT) from vial 1 (SOURCE)	Add 1 μ l of derivatization reagent from vial #1 in the autosampler tray to the derivatization reagent and sample already in the needle capillary.
7	MIX	6 μ l (AMOUNT) in seat (SOURCE), at 500 μ l/min (SPEED), 8 times (REPEAT)	Mix 6 μ l with a repetition of 8 times at a speed of 500 μ l/minute in the seat.
8	WAIT	0.5 minutes (WAIT)	Wait for 0.5 minute for the mixture to completely react.
9	INJECT		Inject the mixture and start the analysis.

A

- action key, 23
 - active signal, 47
 - add button, 35, 48
 - adjusting view, 35
 - Agilent ChemStation, 51
 - features, 58
 - restrictions, 58
 - Agilent on internet, 145
 - align transport arm, 79
 - alphabetic characters, 23
 - alphanumeric characters, 23, 39
 - alphanumeric keypad, 24
 - analog output
 - diode array detector, 92
 - fluorescence light detector, 92
 - variable wavelength detector, 92
 - analysis
 - automated, 44
 - gradient, 109
 - isocratic, 103
 - multiple-vial, 117
 - start, 108, 116
 - time programming, 43
 - analysis screen, 25
 - displayed modules, 35
 - information, 25
 - menu structure, 26
 - printing, 49
 - settings, 26
 - setup view, 35
 - APG Remote
 - interface, 51
 - power on, 51
 - prepare, 52
 - ready, 52
 - shut down, 51
 - start, 52
 - start request, 52
 - stop, 51
 - automated analyses, 44
 - auto-on, 36
 - autosampler
 - align transport arm, 79
 - auxiliary signals, 78
 - configuration, 76
 - defaults, 80
 - draw position offset, 77
 - draw speed, 77
 - eject speed, 77
 - entering settings, 76
 - external contacts, 76
 - injection modes, 77
 - injection volume, 76
 - injector program, 76, 77
 - interfaces, 76
 - missing vial, 76
 - needle seat volume, 76
 - needle wash, 77
 - optimization, 77
 - overlapped injections, 77
 - prefetch, 77
 - resetting, 80
 - runtimes, 76
 - sample throughput, 77
 - shortening analysis time, 77
 - syringe volume, 76
 - tests, 80
 - thermostat, 78
 - theta correction, 79
 - timetable, 76
 - tracking history, 79
 - trays, 76
 - troubleshooting, 80
 - using, 75
 - wash mode, 77
 - wash vial, 77
 - x correction, 79
 - auxiliary signals, 78
 - available modules, 35
 - available signals, 46
- ## B
- backspace, 40
 - balance, 90
 - balance button, 90
 - banner text, 49
 - baseline resetting, 90
 - baudrate, 49
 - BCD interface, 51
 - boot loader firmware, 58
 - buttons, 65
 - add, 35, 48
 - align, 79
 - analog, 93
 - balance, 90
 - calibration, 44
 - column ID, 97
 - column switch, 98
 - configure, 36
 - copy >>, 42
 - cursor, 47
 - delete, 44
 - EMF, 49
 - end actions, 44
 - execute, 55
 - function, 21
 - fw update, 56
 - identify, 48
 - initialize, 42
 - insert, 43, 119
 - interfaces, 51, 76
 - lamp, 88
 - logbook, 44
 - menu, 21
 - method, 26, 38
 - module, 38
 - move, 46
 - on/off, 26, 35, 64, 72, 88, 96
 - page, 49
 - PC card, 42
 - plot, 26
 - previous, 22
 - print, 50
 - printer, 49
 - records, 48
 - related, 21
 - remove/add, 35
 - rescale, 31, 47, 115
 - reset, 69, 80
 - restore, 47
 - right, 97
 - save as, 39
 - search, 21
 - select, 46
 - select file, 55
 - sequence, 26, 44
 - serial, 49
 - settings, 25, 26, 39
 - setup, 35, 46, 50
 - signal, 80
 - spooler, 50
 - start, 126
 - start/stop, 26

temp, 96
tests, 68, 79, 80, 91, 99
timetable, 26, 43
transfer, 55
views, 24, 25, 26
write, 97
zero, 90

C

calibration, 44, 120, 121
 detector, 92
 multiple-level, 121
 sequence, 120, 121
 single-level, 120
 wavelength, 92
calibration button, 44
CAN (Controller Area Network), 34
changing solvents, 73
channels, 65
chromatogram, 114, 115
coexecution, 58
column, 104
column compartment
 column, 104
 column ID, 99
 column identification, 97
 column switching valve, 98
configuration, 97
 heater modes, 98
 ID tag, 97
 modes, 98
 on/off, 96
 resetting, 100
 settings, 96
 tests, 99
 tracking history, 99
 tracking ID history, 99
 troubleshooting, 99
 using, 95
column ID button, 97
column ID module, 97
column switch button, 98
column switching valve, 98
compressibility, 66
configure, 36
context menu, 22
contrast, 35
control module

Agilent ChemStation, 58
available modules, 35
coexecution, 58
contrast, 35
display, 35
displayed modules, 35
features, 19
firmware update, 53
general functions, 35
installing, 34
keys, 20
overview, 20
product number, 20
restrictions, 58
serial number, 20
software, 24
troubleshooting, 36
 user interface, 24
 working with, 33
Control Module Versions, 31
 G1323A, 31
 G1323B, 31
cursor, 47
cursor button, 47

D

date and time, 35
degasser
 changing solvents, 73
 on/off, 72
 removing gas bubbles, 72
 starting, 72
 status lamp, 72
 using, 71
deleting a method, 41
deleting all methods, 41
detector
 baseline resetting, 90
 diode array, 87
 fluorescence light, 87
 variable wavelength, 87
diode array detector, 87
 analog outputs, 90
 autobalance, 88
 balance, 90
 bandwidth, 88
 interfaces, 90
 lamp on/off, 88

peakwidth, 88
resetting, 92
resetting the baseline, 90
settings, 88
slit width, 88
tests, 91
tracking history, 91
troubleshooting, 91
 wavelength, 88
direction keys, 23
display, 46
 contrast, 35
 settings, 35
displayed modules, 35
dot (.) key, 40
draw position offset, 77
draw speed, 77

E

early maintenance feedback (EMF), 19, 28, 68
 counters, 55
 indicator, 48
 logbook, 49
 setting limits, 49
eject speed, 77
EMF button, 49
end actions, 44
end actions button, 44
enter key, 22
error logbook, 48
esc key, 22
execute button, 55
external devices, 51

F

features, 19
firmware
 checking version, 56
 operation mode, 56
 resident mode, 56
 version number, 48, 56
firmware update, 53
 boot loader firmware, 58
 checking version, 56
 control module, 53
 execute, 55
 LC module, 54

-
- operation firmware, 57
 - preparing, 56
 - resident, operation, 54
 - select file, 55
 - switching modes, 56
 - transfer, 55
 - version number, 56
 - fluorescence light detector, 87
 - 3D scan, 89
 - analog outputs, 90
 - emission wavelength, 89
 - excitation wavelength, 89
 - history, 91
 - interfaces, 90
 - lamp on/off, 88
 - multi wavelength, 89
 - peakwidth, 89
 - phosphorescence detection, 89
 - PMT-Gain, 89
 - resetting, 92
 - resetting the baseline, 90
 - settings, 89
 - troubleshooting, 91
 - fluorescence light detectors
 - tests, 91
 - function buttons, 21
 - G**
 - general functions, 35
 - GLP compliance, 19
 - GPIO interface, 51
 - gradient analysis, 109, 111, 115
 - graphical display, 46
 - graphs, 46
 - gripper, 79
 - H**
 - help, 21
 - history
 - autosampler, 79
 - column compartment, 99
 - detector, 91
 - pump, 68
 - I**
 - i (info) key, 21
 - icons
 - EMF, 28, 48
 - ID tag, 97
 - identify button, 48
 - identify modules, 48
 - initialize, 42
 - injection mode, 77
 - injection valve
 - manual, 81
 - injection volume, 76
 - injector program, 77, 129
 - creating, 130
 - entering settings, 130
 - example, 132
 - printing, 50
 - insert button, 119
 - interfaces, 51, 76
 - APG remote, 51
 - BCD, 51
 - GPIO, 51
 - MIO, 51
 - serial/RS232, 51
 - interfaces button, 51
 - internal diagnostics, 36
 - internet, 145
 - isocratic analysis, 103
 - isocratic standard sample, 104
 - K**
 - keypad, 24
 - keys
 - action, 23
 - alphanumeric, 23
 - direction, 23
 - enter, 22
 - esc, 22
 - info, 21
 - keypad, 24
 - menu, 22
 - navigation, 23
 - numeric, 23
 - overview, 20
 - selection, 23
 - L**
 - lamp button, 88
 - lamp ignition routine, 92
 - lamp on/off
 - diode array detector, 88
 - fluorescence light detector, 88
 - variable wavelength detector, 88
 - LC system
 - configuring, 36
 - firmware update, 54
 - general functions, 36
 - shut-down, 44
 - LED, 48
 - liquimeter, 68
 - logbook button, 44
 - logbooks, 48
 - column compartment, 99
 - detector, 91
 - error, 48
 - maintenance, 48
 - print, 36, 50
 - sequence, 44
 - setting up, 48
 - system, 48
 - M**
 - m (menu) key, 22, 40
 - maintenance, 19, 48
 - maintenance logbook, 48
 - manual injection valve, 81
 - filling loops, 83
 - filling sample loop, 83, 85
 - filling sample loop partially, 83
 - making an injection, 82
 - positions, 82
 - sample, 84
 - margins, 49
 - maximize plot screen, 47
 - menu buttons, 21
 - method, 26
 - deleting, 41
 - deleting all, 41
 - exchanging with PC card, 41
 - loading, 38
 - modifying, 39
 - partial method, 38
 - PC card, 38
 - PC card load, 41
 - printing, 50
 - protecting, 40
 - save to PC card, 41
 - saving, 40, 106, 114
 - specifying a name, 39
 - types, 38
-

user ID, 41
 working with, 38
 method button, 38
 MIO interface, 51
 missing vial, 76
 modifying a method, 39
 module button, 38
 Module Status, 26
 move button, 46
 multiple wavelength detector
 autobalance, 89
 bandwidth, 89
 peakwidth, 89
 slit width, 89
 multiple wavelength detector
 settings, 89
 multiple-level calibration, 121
 multiple-level recalibration
 multiple standards, 123
 multiple-vial analysis, 117

N
 navigation key, 23
 needle seat volume, 76
 needle wash, 77
 numeric keypad, 24
 numeric keys, 23

O
 on error, 36
 on/off, 65
 column compartment, 96
 degasser, 72
 diode array detector, 88
 fluorescence light detector, 88
 pump, 64
 variable wavelength detector, 88
 on/off button, 26, 35, 65, 88, 96
 online help, 21
 online information system, 21
 operation firmware, 54
 operation mode, 56
 operation time, 48
 optimizing autosampler, 77
 overlapped injections, 77

P
 page button, 49

page layout, 49
 paper size, 49
 parameter changes, 48
 partial method, 38
 PC card
 initialize, 42
 inserting and removing, 37
 method exchanging, 41
 method loading, 41
 method saving, 41
 PC card button, 42
 piston stroke, 67
 plot button, 26
 plot screen, 30, 46
 active signal, 47
 cursor, 47
 display exact values, 47
 maximize, 47
 print, 50
 print graph, 47
 rescale, 46
 signal unit axis, 47
 time axis, 46
 toggling signals, 47
 x / y values, 47
 x axis, 46
 y axis, 47
 pop-up menu, 24
 power on signal, 51
 prefetch, 77
 prepare signal, 52
 previous button, 22
 print
 delete job, 50
 graph, 47
 interface, 51
 logbooks, 50
 number of copies, 50
 plot, 50
 print button, 50
 printer button, 49
 printer setup, 49
 printing screens, 49
 product number, 20, 48
 protecting a method, 40
 pump, 63
 adjusting compressibility, 66
 channels and composition, 65, 112

 on/off, 64
 piston, 67
 purging, 64
 resetting, 69
 resetting settings, 69
 settings, 64
 shut-down, 67
 stroke, 67
 tracking history, 68
 troubleshooting, 68
 using, 63
 valve, 64
 purge
 procedure, 65
 pump, 64
 valve, 64

R
 ready signal, 52
 recalibration, 121
 alter/multi, 44
 recalibration with same group of stan-
 dards, 121
 records nutton, 48
 refractive index detector
 automatic recycling, 89
 cell temperature, 89
 peakwidth, 89
 polarity, 89
 settings, 89
 zero, 90
 related button, 21
 remove button, 35
 rescale button, 31, 47, 115
 rescale plot screen, 46
 reset button, 69, 80
 resetting
 autosampler, 80
 autosampler defaults, 80
 baseline, 90
 column compartment, 100
 detector, 92
 pump, 69
 pump settings, 69
 resident firmware, 54
 resident mode, 56
 restore button, 47
 right button, 97

RS232 interface, 51
running an analysis, 103
running injector program, 129

S

sample
 amount, 84
sample loop
 completely filling, 83
samples screen, 30
 printing, 49
save as button, 39
saving a method, 40
saving method, 106, 114
screens
 analysis, 24, 25, 26
 config, 49
 method, 50
 plot, 24, 30, 46
 program, 50
 records, 48
 samples, 24, 30
 sequence, 44
 status, 24, 29
 system, 24, 28
 timetable, 39
 views, 26
search button, 21
select button, 46
select file button, 55
select signals, 46
selection keys, 23
sequence, 26, 44, 119, 120, 121
 calibration, 44, 120, 121
 end actions, 44
 logbook, 44
 printing, 50
 recalibration, 44
sequence button, 44
sequence modes, 126
serial button, 49
serial interface, 51
serial number, 20, 48
settings
 detectors, 88
 diode array detector, 88
 fluorescence light detector, 89
 multiple wavelength detector, 89

 refractive index detector, 89
 settings button, 25, 39
 setup button, 35, 46, 50
 setup logbooks, 48
 setup view, 35
 shut down signal, 51
 shut-down pump, 67
 signal unit axis, 47
 signals button, 80
 single-level calibration, 120
 solvent composition, 65, 104
 solvents, 73
 spooler button, 50
 standard isocratic sample, 104
 standard mode, 126
 start button, 126
 start request signal, 52, 126
 external, 127
 start signal, 52, 126
 start/stop button, 26
 starting an analysis, 108, 116
 status screen, 29
 stop signal, 51
 stroke volume, 67
 syringe volume, 76
 system logbook, 48
 system screen
 information, 28
 menu structure, 29

T

tag, 97
tests
 autosampler, 79, 80
 column compartment, 99
 detector, 91
tests button, 68, 79, 80, 99
thermostatted autosampler, 78
 ambient temperature, 78
 auxiliary signals, 78
 display signals, 78
 heat sink, 78
 heat sink fan, 78
 parameters, 78
 plot, 78
 settings, 78
 signals, 78
theta correction, 79

time and date, 35
time axis, 46
time programming, 43
timeout, 36
timetable, 26, 43
 autosampler, 76
 copy / paste, 44
 printing, 50
 setting, 43
timetable button, 43
toggling signals, 47
transfer, 56
transfer button, 55
transport arm aligning, 79
trays, 76
troubleshooting
 autosampler, 80
 column compartment, 99
 control module, 36
 detectors, 91
 logbook, 36
 pump, 68
turning on/off
 column compartment, 96

U

units, 36
updating firmware, 53
user ID, 41

V

valve
 column compartment, 98
 filling sample loop, 83, 85
 filling sample loop partially, 83
 making an injection, 82
 manual, 81
 opening, 64
 positions, 82
 sample, 84
variable wavelength detector, 87
 analog output, 90
 balance, 90
 interfaces, 90
 lamp ignition routine, 92
 lamp on/off, 88, 92
 resetting, 92
 resetting the baseline, 90

- settings, 88
- signal polarity, 88
- spectrum, 88
- tests, 91
- tracking history, 91
- troubleshooting, 91
- wavelength, 88
- wavelength calibration, 92
- version number, 56
 - checking, 56
- views, 26
- views button, 24, 25

W

- warranty
 - responsibility of Agilent Technologies, 141
 - services, 141
 - statement, 139
- wash mode, 77
- wash vial, 77
- washing the needle, 77
- wavelength calibration, 92
- wear count, 68
- working with control module, 33
- working with methods, 38
- write button, 97

X

- x / y values, 47
- x axis, 46
- x correction, 79

Y

- y axis, 47

Z

- zero button, 90

Warranty Statement

All Chemical Analysis Products

Agilent Technologies warrants its chemical analysis products against defects in materials and workmanship. For details of the warranty period in your country, call Agilent. During the warranty period, Agilent will, at its option, repair or replace products which prove to be defective. Products that are installed by Agilent are warranted from the installation date, all others from the ship date.

If buyer schedules or delays installation more than 30 days after delivery, then warranty period starts on 31st day from date of shipment (60 and 61 days, respectively for products shipped internationally).

Agilent warrants that its software and firmware designed by Agilent for use with a CPU will execute its programming instructions when properly installed on that CPU. Agilent does not warrant that the operation of the CPU, or software, or firmware will be uninterrupted or error-free.

Limitation of Warranty

Onsite warranty services are provided at the initial installation point. Installation and onsite warranty services are available only in Agilent service travel areas, and only in the country of initial purchase unless buyer pays Agilent international prices for the product and services. Warranties requiring return to Agilent are not limited to the country of purchase.

For installation and warranty services outside of Agilent's service travel area, Agilent will provide a quotation for the applicable additional services.

If products eligible for installation and onsite warranty services are moved from the initial installation point, the warranty will remain in effect only if the customer purchases additional inspection or installation services, at the new site.

The foregoing warranty shall not apply to defects resulting from:

- 1 improper or inadequate maintenance, adjustment, calibration, or operation by buyer,
- 2 buyer-supplied software, hardware, interfacing or consumables,
- 3 unauthorized modification or misuse,
- 4 operation outside of the environmental and electrical specifications for the

Warranty Statement

product,

- 5 improper site preparation and maintenance, or
- 6 customer induced contamination or leaks.

THE WARRANTY SET FORTH IS EXCLUSIVE AND NO OTHER WARRANTY, WHETHER WRITTEN OR ORAL, IS EXPRESSED OR IMPLIED. AGILENT SPECIFICALLY DISCLAIMS THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE.

Limitation of Remedies and Liability

THE REMEDIES PROVIDED HEREIN ARE BUYER'S SOLE AND EXCLUSIVE REMEDIES. IN NO EVENT SHALL AGILENT BE LIABLE FOR DIRECT, INDIRECT, SPECIAL, INCIDENTAL, OR CONSEQUENTIAL DAMAGES (INCLUDING LOSS OF PROFITS) WHETHER BASED ON CONTRACT, TORT OR ANY OTHER LEGAL THEORY.

Responsibilities of the Customer

The customer shall provide:

- 1 access to the products during the specified periods of coverage to perform maintenance,
- 2 adequate working space around the products for servicing by Agilent personnel,
- 3 access to and use of all information and facilities determined necessary by Agilent to service and/or maintain the products (insofar as these items may contain proprietary or classified information, the customer shall assume full responsibility for safeguarding and protection from wrongful use),
- 4 routine operator maintenance and cleaning as specified in the Agilent operating and service manuals, and
- 5 consumables such as paper, disks, magnetic tapes, ribbons, inks, pens, gases, solvents, columns, syringes, lamps, septa, needles, filters, frits, fuses, seals, detector flow cell windows, and so on.

Responsibilities of Agilent Technologies

Agilent Technologies will provide warranty services as described in Table 5.

Table 5

Warranty Services

Services During Warranty*	Warranty Period**	Type
Agilent 1100 Series of Modules	1 Year	Onsite
GC, LC, UV-Visible, and LAS supplies and accessories	90 Days	Onsite
Columns and Consumables***	90 Days	Return to Agilent
Gas Discharge and Tungsten Lamps	30 Days	Return to Agilent
Repairs performed on-site by Agilent****	90 Days	Onsite

* This warranty may be modified in accordance with the law of your country. Please consult your local Agilent office for the period of the warranty, for shipping instructions and for the applicable wording of the local warranty.

** Warranty services are included as specified for Analytical products and options purchased concurrently provided customer is located within a Agilent defined travel area. Agilent warranty service provides for 8 a.m. to 5 p.m. on-site coverage Monday through Friday, exclusive of Agilent holidays.

*** Columns and Consumables are warranted to be free from defects for a period of 90 days after shipment and will be replaced on a return-to-Agilent basis if unused.

**** Agilent repair warranty is limited to only the item repaired or replaced.

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.

General

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 fuseholders 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 as much as possible. When inevitable, this should 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.

Safety Information

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.

Safety Symbols

Table 6 shows safety symbols used on the instrument and in the manuals.

Table 6

Safety Symbols	
Symbol	Description
	The apparatus is marked with this symbol when the user should refer to the instruction manual in order to protect the apparatus against damage.
	Indicates dangerous voltages.
	Indicates a protected ground terminal.
	Eye damage may result from directly viewing the light produced by the Xenon flash lamp used in this product. Always turn the xenon flash lamp off before removing it.

WARNING

A warning alerts you to situations that could cause physical injury or damage to the equipment. Do not proceed beyond a warning until you have fully understood and met the indicated conditions.

Radio Interference

CAUTION

A caution alerts you to situations that could cause a possible loss of data. Do not proceed beyond a caution until you have fully understood and met the indicated conditions.

Radio Interference

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

Test and Measurement

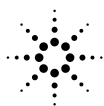
If test and measurement equipment is operated with equipment unscreened cables and/or used for measurements on open set-ups, the user has to assure that under operating conditions the radio interference limits are still met within the premises.

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”



Agilent Technologies

Innovating the HP Way

In This Book

This book provides information about the control module and how to operate the Agilent 1100 Series modules and system for HPLC.

- The Agilent 1100 Series Control Module
- Working with the Control Module
- Using the Pump
- Using the Degasser
- Using the Autosampler
- Using the Manual Injection Valve
- Using the Detectors
- Using the Column Compartment
- Running an Isocratic Analysis
- Running a Gradient Analysis
- Running Multiple-Vial Analyses
- Running an Injector Program



G1323-90006