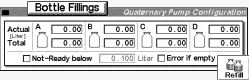
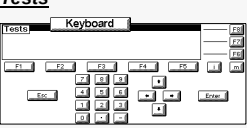
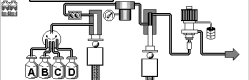
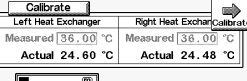
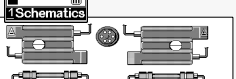
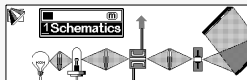
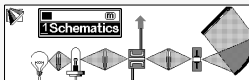
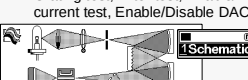


LC System	Quat Pump	Autosampler	Column Compartment	Diode Array Detector	Multiple Wavelength Detector	Variable Wavelength Detector
Settings Runtimes Stoptime [min] Posttime [min] <i>Module runtime settings override LC settings</i>      	Settings Solvent Composition %B, %C, %D $%A = 100 - \%B - \%C - \%D$ Flow [ml/min] More... Flow Ramp [ml/min ²] Compressibility [10 ⁻⁹ /bar] Primary Channel [A,B,C,D,Auto] Stroke [μl] Pressure Lower/Upper Pressure Limits [bar]         	Settings Injection Volume [μl] ☒ Standard Injection ☐ with needle wash ☒ Injector Program Default wash vial More... Settings Draw Speed [μl/min] Eject Speed [μl/min] Draw Position Offset [mm] More... Optimization Prefetch Vial / Overlap Injection Cycle at elapsed runtime [min] ☒ always keep transport arm next to last vial        	Settings Temperature Left [°C] Temperature Right [°C] ☒ enable left/right separate mode More... Enable Analysis ☒ with any temperature ☒ temp. within setpoint +/- [°C]         	Settings Sample/Ref. Wavelengths (A, B) [nm] Sample/Ref. Bandwidths (A, B) [nm] Peakwidth(Responsetime) [min] Slit [nm] More... Settings Auto Balance: ☒ Prerun/ ☒ Postrun Margin for negative Absorbance [mAU] Required Lamps ☒ UV/ ☒ VIS More... Signals C-E Sample/Ref. WL & BW (C, D, E) [nm] Analogs Out 1/2: Zero Offset [%] Out 1/2: Attenuation [mAU] Spectrum Range [nm] Store Spectra Threshold [mAU]          	Settings Sample/Ref. Wavelengths (A, B) [nm] Sample/Ref. Bandwidths (A, B) [nm] Peakwidth(Responsetime) [min] Slit [nm] More... Settings Auto Balance: ☒ Prerun/ ☒ Postrun Margin for negative Absorbance [mAU] Required Lamps ☒ UV/ ☒ VIS More... Signals Sample/Ref. Bandwidths (C, D, E) [nm] Analogs Out 1/2: Zero Offset [%] Out 1/2: Attenuation [mAU] Spectrum Range [nm]         	Settings Wavelength [nm] Peakwidth (Responsetime) [min] Signal Polarity More... Auto Balance: ☒ Prerun/ ☒ Postrun Margin for negative Absorbance [mAU] ☒ Not-Ready when lamp is off Analogs Signal Source Zero Offset [%] Attenuation [mAU] Spectrum Range[nm]    
Control    	Control    	Control       	Control    	Control     	Control     	Control  
Configure Auto-On [month/day/time] ☒ turn on Module/LC System ☒ load Method After Error Condition: ☒ turn off Module/LC System ☒ run Method      	Configure   	Configure Syringe Volume [μl] Seat Capillary [μl] Multiple Draw Wait [s] On Missing Vial  	Configure ☒ Disable leak detection ☒ At Power-On turn temperature on Column ID Stationary Phase Geometry; Particle Size; Void Volume Product -, Serial -, Batch Number Write Column ID More... Max Pressure, - Temperature, - pH Number of Injections Comment   	Configure Options Analog Voltage Range Out 1 Analog Voltage Range Out 2 At Power-On turn ☒ UV/ ☒ VIS lamp on  	Configure Options Analog voltage range ☒ At Power-On turn UV lamp on Lamp Type  	Configure Options Analog voltage range ☒ At Power-On turn UV lamp on Lamp Type  
Tests    	Tests         	Tests Functions: Valve bypass/mainpass, Syringe home/draw, Needle up into vial, Fetch/Return vial; Align Tray Tray alignment/Vial teaching Align Transport X Correction [mm] Theta Correction [deg] Change Change needle/piston/arm          	Tests         	Tests Functions: Start/Stop test chromatogram, Enable/Disable DAC test pattern            	Tests Functions: Start/Stop test chromatogram, Enable/Disable DAC test pattern            	Tests Functions: Start/Stop test chromatogram, Grating test, Filter test, Enable/Disable Dark current test, Enable/Disable DAC test pattern          

For information on other Agilent 1100 modules, please refer to the Modules Reference Manual.

Samples Screen

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

☒ **Vial Range** Vial 1 to 8 #Inj 1

Method TEST Stoptime 5.00 min

☐ **Sequence** Line 1 to 8

End Actions Timetable Method Sequence Views

On end of ...
☐ Load Method *ACTUAL*
☒ Turn off Nothing
☐ Turn off complete HPLC system
 Timeout 0.5 min

1 Balance
2 Print Screen
3 Restart

Start On/Off Plot

Status Screen

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

Flow 0.300 ml/min Temp 25.80 °C

% B 20.0 % 1.117 mA

12 bar | Ripple -0.3 % λ 250 nm

Elapsed 0.00 min

Start On/Off Plot

1 Balance

System Screen

System Pump ☒ Temp ☒ Lamp ☒ Time 0.00 Idle Ready *

Fri 15:20

1 Date & Time
2 Save Logbook
3 Print Logbook
4 Setup Logbook
5 Restart

Module	Message	Id	Date	Time
Quat Pump	Setpoint changed	INF	03/26	15:19:26
Quat Pump	Pump on	STAT	03/26	15:18:46
VW Detector	UV lamp burn time limit	EMF	03/26	15:18:13
Col Comp	Calibration done	STAT	03/26	15:18:09
MW Detector	VIS lamp on	STAT	03/26	15:18:13

Control Configure Tests Records Views

Start On/Off Plot



Agilent Technologies
Innovating the HP Way



G1323-90007

© Copyright Agilent Technologies 2000
Printed in Germany 01/00
Part No. G1323-90007

Analysis Screen

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

Fri 15:21 1 - 1 # 1 TEST 5.00

Flow 0.300 ml/min 0.3 µl L OFF °C A λ BW 1 0 On/Off Plot

0.300 0.300 25.81 °C Samp 250/100

16 bar 0.0 µl 24.73 °C Ref 360/100 23.892 °C

Settings Timetable Method Sequence Views

1 Pump On/Off
2 Bottle Fillings
3 Reset
4 Status
5 Setup View
6 Restart

1 Program
2 Reset
3 Status
4 Setup View
5 Restart

1 Temp On/Off
2 Column ID
3 Reset
4 Status
5 Setup View
6 Restart

1 Lamp On/Off
2 Balance
3 Reset
4 Status
5 Setup View
6 Restart

1 Date & Time
2 Print Screen
3 About
4 Setup View
5 Restart

Start On/Off Plot

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

Module Setting Value

Quat Pump Autosampler Col Comp VW Detector DIA Detector FL Detector

Information stored in Methods:
 - Runtimes - User Information
 - Settings - Injector Program
 - Timetable - Optimization

Save As Default Print

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

Line Vials # Inj.Vol Method Wait Calibration

Information stored in Sequences:
 - Vial Range Information
 - Injection Volume - Wait Time (after Method Load)
 - Method Name - (Re-) Calibration Settings

Start Insert Delete

End Actions PC-Card Method Logbook

