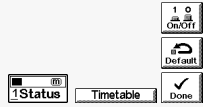
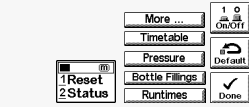
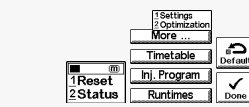
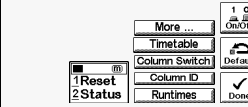
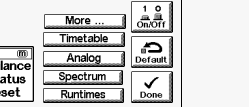
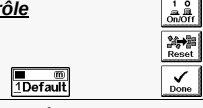
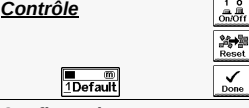
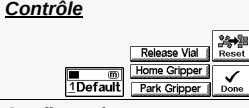
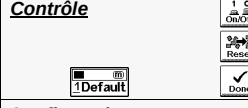
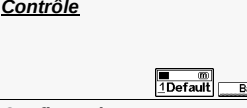
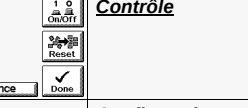
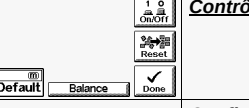
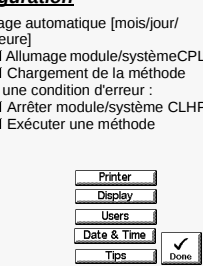
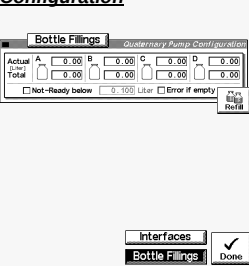
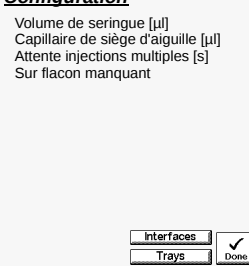
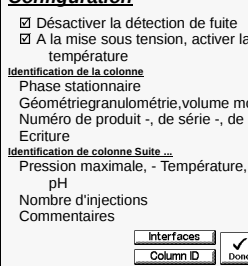
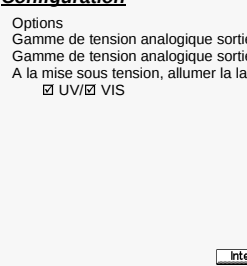
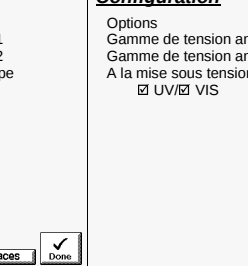
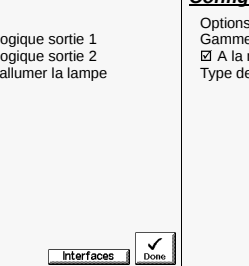
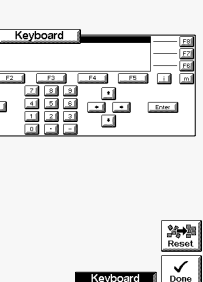
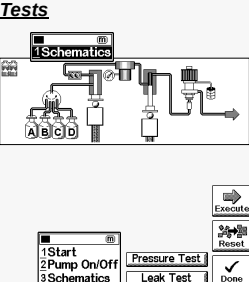
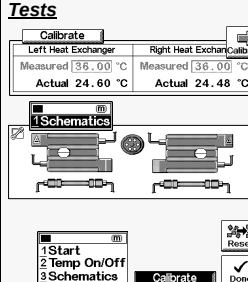
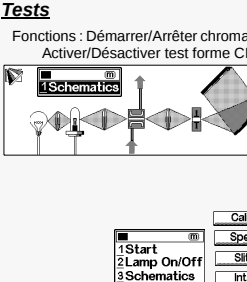
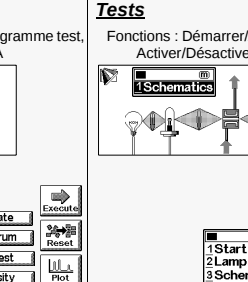
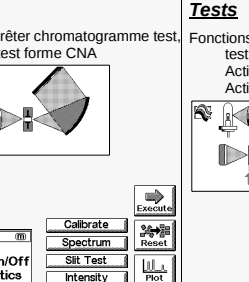


Système CLHP	Pompe quaternaire	Passeur automatique d'échantillons	Compartment à colonne	Détecteur à barrette de diodes	Détecteur multi-longueur d'onde	Détecteur à longueur d'onde variable
Paramétrages (settings) Runtimes (temps d'analyse) Stoptime [min] Temps avant arrêtPosttime [min] isotherme post-analytique Paramétrages des temps d'analyse du module Forçage des paramétrages CLHP 	Paramétrages (settings) Solvent Composition (composition du solvant) %B, %C, %D $%A = 100 - \%B - \%C - \%D$ Flow [ml/min] Débit [ml/min] Suite ... Flow Ramp - rampe de débit [ml/min ²] Compressibility - compressibilité [10 ⁻⁹ /bar] Primary Channel - voie primaire [A,B,C,D,Auto] Stroke - course du piston [μl] Pressure - pression Lower/Upper Pressure Limits - Limites supérieure/inférieure de pression [bar] 	Paramétrages (settings) Volume d'injection [μl] ☒ Injection normale ☒ avec rinçage d'aiguille ☒ Programme d'injecteur rinçage de flacon par défaut Suite ... Paramétrages Vitesse de prélèvement [μl/min] Vitesse d'expulsion [μl/min] Décalage de position de prélèvement [mm] Suite ... Optimisation Pré-sélection du flacon/Superposition cycle d'injection au temps d'analyse écoulé [min] ☒ toujours conserver le bras près du dernier flacon 	Paramétrages (settings) Température Gauche [°C] Injection normale Température Droit [°C] ☒ activer mode distinct gauche/droite Suite ... ☒ Activer Analyse ☒ à n'importe quelle température ☒ temp. dans la limite du point de consigne +/- [°C] 	Paramétrages (settings) Echantillon/Ref. Longueurs d'onde (A, B) [nm] Echantillon/Ref. Bande passantes (A, B) [nm] Largeur de pic(Responsetime) [min] Fente [nm] Suite ... Paramétrages Stabilisation : ☒ pré-analytique/ ☒ post-analyse Marge de l'absorbance négative [mDO] Lampes nécessaires ☒ UV/ ☒ VIS Suite ... Signaux C-E Echantillon/Ref. Long. onde & Bande passante (C, D, E) [nm] Analog Sortie 1/2 : décalage zéro [%] Sortie 1/2 : Atténuation [mDO] Spectre Gamme [nm] Stockage des spectres Seuil [mDO] 	Paramétrages (settings) Echantillon/Ref. Longueurs d'onde (A, B) [nm] Echantillon/Ref. Bandes passantes (A, B) [nm] Largeur de pic(Responsetime) [min] Fente [nm] Suite ... Paramétrages Stabilisation : ☒ pré-analytique/ ☒ post-analyse Marge de l'absorbance négative [mDO] Lampes nécessaires ☒ UV/ ☒ VIS Suite ... Signaux Echantillon/Ref. Bandes passantes (C, D, E) [nm] Analog Sortie 1/2 : décalage zéro [%] Sortie 1/2 : Atténuation [mDO] Spectre Gamme [nm] 	Paramétrages (settings) Longueur d'onde [nm] Largeur de pic (Responsetime) [min] Polarité du signal Suite ... Stabilisation : ☒ pré-analytique/ ☒ post-analyse Marge de l'absorbance négative [mDO] ☒ Non-Prêt lorsque la lampe est éteinte Analog Source du signal Décalage zéro [%] Atténuation [mDO] Spectre Gamme [nm] 
Contrôle 	Contrôle 	Contrôle 	Contrôle 	Contrôle 	Contrôle 	Contrôle 
Configuration Allumage automatique [mois/jour/heure] ☒ Allumage module/système CPL ☒ Chargement de la méthode Après une condition d'erreur : ☒ Arrêter module/système CLHP ☒ Exécuter une méthode 	Configuration 	Configuration Volume de seringue [μl] Capillaire de siège d'aiguille [μl] Attente injections multiples [s] Sur flacon manquant 	Configuration ☒ Désactiver la détection de fuite ☒ A la mise sous tension, activer la température Identification de la colonne Phase stationnaire Géométriegranulométrie, volume mort Numéro de produit -, de série -, de lot Ecriture Identification de colonne Suite ... Pression maximale, - Température, - pH Nombre d'injections Commentaires 	Configuration Options Gamme de tension analogique sortie 1 Gamme de tension analogique sortie 2 A la mise sous tension, allumer la lampe ☒ UV/ ☒ VIS 	Configuration Options Gamme de tension analogique sortie 1 Gamme de tension analogique sortie 2 A la mise sous tension, allumer la lampe ☒ UV/ ☒ VIS 	Configuration Options Gamme de tension analogique ☒ A la mise sous tension, allumer la lampe UV Type de lampe 
Tests 	Tests 	Tests Fonctions : vanne contournement/circuit principal, Seringue position de départ/prélèvement, Aiguille en l'air/dans le flacon, Flacon : prendre/Remettre en place ; Alignement du plateau Alignement du plateau/reconnaissance de flacon Alignement du passeur Correction X [mm] Correction Theta [deg] Changement Changement aiguille/piston/bras 	Tests 	Tests Fonctions : Démarrer/Arrêter chromatogramme test, Activer/Désactiver test forme CNA 	Tests Fonctions : Démarrer/Arrêter chromatogramme test, Activer/Désactiver test forme CNA 	Tests Fonctions : Démarrer/Arrêter chromatogramme test, test du réseau holographique, test de filtre, Activer/Désactiver test de courant d'obscurité, Activer/Désactiver test forme CNA 

Pour plus d'information sur les autres modules de la Série Agilent 1100, veuillez consulter les manuels de références des différents modules.

Ecran Echantillons

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 ...
 Vial Range
 Timeout 0.5 min

Load Method ACTUAL *
 Turn off Nothing
 Turn off complete HPLC system

1 Balance
 2 Print Screen
 3 Restart

Start
 On/Off
 Plot

Ecran Etats

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

Ecran Système

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

Fri 15:20

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

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

Start
 On/Off
 Plot



Agilent Technologies



G1323-93007

© Copyright Agilent Technologies 2000
 Imprimé en Allemagne 01/00
 Réf. G1323-93007

Ecran Analysis

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

25.81 °C Samp 250/100

R 20.00 °C Ref 360/100

16 bar 0.0 µl 24.73 °C 23.892 °C

Settings Timetable Method Sequence Views

1 Pump On/Off
 2 Bottle Filling
 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	Information stockées dans méthodes
Autosampler	- Temps d'analyse Information utilisat
Col Comp	- Paramétrages - Program injecteur
VW Detector	- Chronoprogram - Optimisation
DIA Detector	
FL Detector	

Module PC-Card Timetable

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

Line Vials # Inj.Vol Method Wait Calibration

Information stockées dans les séquences:

- Information gamme flacon
- Volume d'injection
- Nom de méthode
- Temps attente (après charg.méthode)
- Paramètres ré-étalonnage

End Actions PC-Card Method Logbook

Start
 On/Off
 Plot

