



708-DS
Dissolution Apparatus
Operator's Manual

Notices

Manual Part Number

70-9058 Rev H
February 2020

Copyright

© Agilent Technologies, Inc. 2020

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.
3501 Stevens Creek Blvd.
Santa Clara, CA 95052 USA

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Content

1	Safety	7
	Electrical Hazards	8
	Warning	9
	Caution	9
	Note	9
	Information Symbols	10
2	Introduction	11
	Conventions Used in this Manual	13
	Serial Number Format	14
3	Setting Up the 708-DS	15
	Initial Setup	16
	Unpacking Procedure	16
	Environmental Requirements for Installation	17
	Clearance	17
	Equipment, Parts, and Accessories As Shipped	18
	Parts and Accessories	18
	Tubing / Cable Connections	19
	708-DS Heater / Circulator Setup	19
	Instrument Level / Water Bath Setup	20
	Level Instrument	20
	Bath Temperature Probe Installation	22
	Handheld Probe Installation	23
	Power Cord Connections	23
	708-DS Initial Power Up	24
	Filling the Water Bath	24
	Level Instrument - Re-verification	26
	Heater / Circulator Power-up	26
	Agilent 708-DS Setup	27
	Vessel Installation	27

Content

Sample Manifold Installation	27
Accessory Installation	31
Setting Basket (Apparatus 1) Heights	37
Setting Paddle (Apparatus 2) Heights	38
4 Operating the 708-DS	39
Main Screen - Options	40
Main Screen - Display Parameters	40
Main Screen - Sample Temperatures	41
Main Screen - Run	42
Main Screen - Stop	48
Main Screen - Drive Unit Up / Down	51
Manual Sampling	51
Main Screen - Manifold Up / Down	53
Main Screen - Lock	53
Main Screen - Unlock or Log In	54
Main Screen - Remote Control	56
Main Screen - Alarms	57
Main Screen - Maintenance Due Icon	57
Menu Screen - System Menu	58
Menu Screen - Method Editor	59
Menu Screen - Reports	67
Menu Screen - Alarms	70
Menu Screen - Instrument	71
Menu Screen - Calibration	76
Menu Screen - Diagnostics	79
Menu Screen - User Access	81
Menu Screen - Mfg. Config	85
General 708-DS Conventions	86
Alphanumeric Data Entry	86
Numeric and Time / Date Entry	87
Parameter Limits / Ranges	88

Content

Agilent 708-DS Notifications	90
Main Screen - Error Conditions	90
Dialog Screens (Green / Blue)	91
Warning Screens (Yellow)	93
Error Screens (Red)	97
5 Maintenance and Troubleshooting	101
Preventive Maintenance	102
Weekly Maintenance	102
Monthly Maintenance	102
Bi-Annual Maintenance	102
Sample Line / Paddle / Basket Care	103
Shafts	103
Baskets	104
Rinsing Sample Lines	105
Rinse Cups (Optional)	106
Vessels and Evaporation Covers	107
Water Bath and Temperature Probes	107
Water Bath Care	107
Water Bath Temperature Probe Accuracy Test	108
Removing or Adjusting Cannulas / Probes	110
Manifold Maintenance	110
Removing the Manifold	110
Accessing the Top Cover	111
Removing Top Cover	111
Replacing the Top Cover	111
Printer Maintenance	112
Removing a Thermal Paper Roll	112
Inserting a Thermal Paper Roll	113
Thermal Printer Test	115
Obtaining Warranty and Other Services	116

Content

This page was intentionally left blank, except for this message.

1

Safety

Electrical Hazards 8

The Agilent 708-DS has been designed and tested so that when used properly you have an accurate, fast, flexible, and safe instrument.

If the equipment is used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

Operation of an Agilent 708-DS involves the use of aqueous liquids and various pharmaceutical dosage forms. Unskilled, improper, or careless use of this instrument can create shock hazards, fire hazards, or other hazards which can cause death, serious injury to personnel, or severe damage to equipment and property.

Information on safety practices is provided with your instrument and operation manuals. Before using your instrument or accessories, you must thoroughly read these safety practices.

Observe all relevant safety practices at all times.

Electrical Hazards

The dissolution apparatus contains electrical circuits, devices, and components operating at dangerous voltages. Contact with these circuits, devices, and components can cause death, serious injury, or painful electric shock.

Panels or covers that are retained by fasteners which require the use of a tool for removal may be opened only by Agilent-trained, Agilent-qualified, or Agilent-authorized service engineers. Consult the manuals or product labels supplied with the dissolution apparatus to determine which parts are operator-accessible.

Application of the wrong supply voltage, connection of the instrument to an incorrectly wired supply outlet, or lack of proper electrical grounding can create a fire hazard or a potentially serious shock hazard and could seriously damage the instrument and any attached ancillary equipment.

Always use a three-wire outlet with ground connection which is adequately rated for the load. The installation must comply with local, state, and federal safety regulations.

Do not connect the instrument to the main power supply until you have made sure that the operating voltage is correctly set for the main power supply in the specific outlet in your laboratory to which the equipment will be connected.

Warning

WARNING

A 'Warning' message appears in the manual when failure to observe instructions or precautions could result in death or injury.

Read all warnings and cautions carefully and observe them at all times.

Caution

CAUTION

A 'Caution' message appears in the manual when failure to observe instructions could result in damage to equipment (Agilent supplied and / or other associated equipment).

Note

NOTE

A 'Note' appears in the manual to give advice or information.

Information Symbols

I

Switches main power on

0

Switches main power off



Indicates single-phase alternating current



Indicates the product complies with the requirements of one or more European Union (EU) directives.



Indicates specific equipment meets standards of safety. These products are safe for use in the workplace for North America.



Indicates that this product must not be disposed of as unsorted municipal waste.

All Agilent products that are subject to the WEEE directive shipped after August 13, 2005 are compliant with the WEEE marking requirements. Such products are marked with the “crossed out wheeled bin” WEEE symbol in accordance with European Standard EN 50419.

For more information on collection, reuse, and recycling systems, please contact your local/regional waste administration, your local distributor, or Agilent.



Indicates the product complies with regulatory compliance requirements of New Zealand and Australia.



2

Introduction

Conventions Used in this Manual 13

Serial Number Format 14

Agilent's 708-DS Dissolution Apparatus is designed for dissolution testing of a variety of pharmaceutical products, including tablets, capsules, and transdermal patches. Interaction with the instrument is simplified with an intuitive, color touchscreen interface. Temperature control is achieved using traditional water bath vessel heating. This apparatus can be tailored to support several innovative features designed to allow unattended dissolution testing, including Dosage Delivery Module (DDM), AutoTemp, and Auto Sampling.

WARNING

The dissolution apparatus contains electrical circuits, devices, and components operating at dangerous voltages. Contact with these circuits, devices and components can cause death, serious injury, or painful electric shock.

CAUTION

Panels or covers that are retained by fasteners which require the use of a tool for removal may be opened only by Agilent-trained, Agilent-qualified, or Agilent-authorized service engineers

Conventions Used in this Manual

- Items you are asked to press are in bold. For example, “press **H** on the keypad”.
- Key sequences you are asked to press appear like this: **MENU > 7**.

Serial Number Format

The serial number contains 10 characters and follows this syntax:

CC1234XXXX

Syntax Code	Meaning	Description
CC	Country of origin	2 alpha characters matching the required trade designation for the country of origin
12	Year of manufacture	'09' for 2009, '10' for 2010, etc.
34	Week of manufacture	'01' for week 1 to '52' for week 52



3

Setting Up the 708-DS

Initial Setup 16

Equipment, Parts, and Accessories As Shipped 18

Tubing / Cable Connections 19

Instrument Level / Water Bath Setup 20

Agilent 708-DS Setup 27

Initial Setup

Complete the following sections to unpack the Agilent 708-DS and ensure it is set up in an appropriate location.

Unpacking Procedure

- 1 Remove the Agilent 708-DS and all other system components from the packing material.

CAUTION

Do not remove the protective foam from the drive unit until it has been raised.

- 2 Inspect the equipment and accessories to ensure there has been no damage during shipment.

Environmental Requirements for Installation

- Humidity: max relative humidity 80% for temperatures up to 31 °C decreasing linearly to 50% relative humidity at 40°C
- Indoor use only
- Pollution Degree: 2
- Installation Category: II
- Altitude: 2000m
- Temperature: 5 °C to 40 °C
- Power: 115/230 V ac, 50/60 Hz, 2.5 A

Main supply voltage fluctuations are not to exceed $\pm 10\%$ of the nominal supply voltage

Clearance

- 1 Prepare the area where the equipment is to be located.
- 2 Ensure a minimum clearance of 30 cm above the unit (when fully raised) and 10 cm at the rear and on both sides of the Agilent 708-DS. Approximately 82 cm x 82 cm total bench space is required per apparatus.
- 3 Orient the Agilent 708-DS and all other system components appropriately.

Equipment, Parts, and Accessories As Shipped

Parts and Accessories

Locate the following items, as applicable, for your system configuration:

708-DS Dissolution Apparatus	Alignment Posts or Evaporation Plugs
Heater / Circulator	Evaporation Covers
6-pin Cable for Heater / Circulator	Blank Vessel Position Covers
Bath Temperature Probe	17-mm Open-end Wrench
External Vessel Temperature Probe	90 ° Allen Key (for Level Adjustment)
Power Cord(s)	Shaft Locking Collars
Manifold Arm Labels	Vessel Centering Tool
6-pin Cable for Heater / Circulator	Printer Paper
Bath Temperature Probe	7/64" T-handle Allen Wrench
Rinse Tray	Stainless Steel Tubing Clamps
Dissolution Vessels	Plastic Tubing - 1/2" outer diameter
Dosage Delivery Modules (DDMs)	25-mm Height Spheres
Basket Shaft Assemblies	Clip-on Basket Height Gauge
Paddle Shaft Assemblies	Bubble Level
Technical Documentation CD	

Tubing / Cable Connections

Complete the following sections to connect the necessary tubing and cables for the Agilent 708-DS.

708-DS Heater / Circulator Setup

CAUTION

Do not turn on the heater / circulator before filling the water bath to avoid damaging the internal heating elements of the circulator.

NOTE

The Agilent Dissolution Heater / Circulator is pre-installed on the apparatus at the factory. If reinstallation or replacement is required, please refer to the *Agilent Dissolution Heater / Circulator Operator's Manual* (located on the Technical Documentation CD) for complete installation instructions.

- 1 Connect the six-pin cable to the rear of the heater / circulator and the other end into the position marked BATH HEATER on the dissolution apparatus rear panel.



Figure 1. BATH HEATER Connector

Instrument Level / Water Bath Setup

Complete the following sections to properly level the instrument and fill the water bath.

Level Instrument

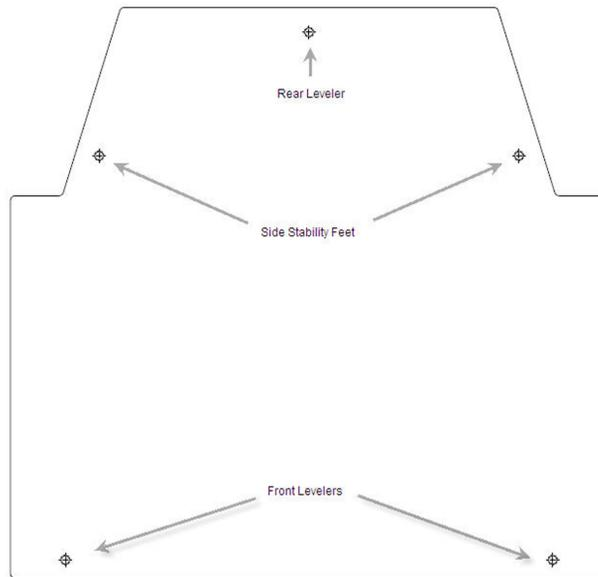


Figure 2. Levelers and Stability Feet

- 1 Remove the black caps from the five (5) screws located on the front, side, and rear of the base plate of the dissolution apparatus.



Figure 3. Level-Adjustment Screw Caps

- 2 Raise the side stability feet located toward the back on either side of the base plate so they are not touching the laboratory bench.

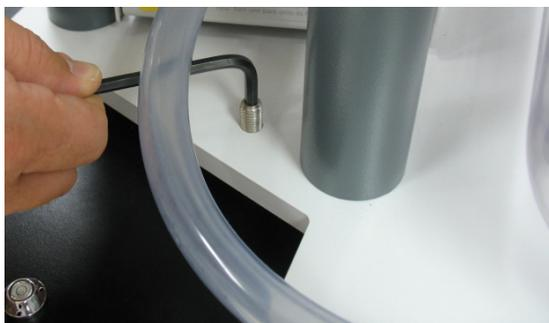


Figure 4. Raising the Stability Feet

- 3 Using the 90° Allen key in the top of the front level-adjustment screws, adjust the screws to achieve left-to-right level within tolerance. It may be necessary to loosen the nut(s) beneath the base plate to allow for adjustment.

- 4 Using the 90° Allen key in the top of the rear level-adjustment screw, adjust the screw to achieve front-to-back level within tolerance. It may be necessary to lift the heater / circulator to gain access to this screw.
- 5 Verify the level of the dissolution apparatus with a bubble or digital level.
- 6 Once the unit is leveled, use the open-end wrench to tighten the nut(s) below the base plate.
- 7 Lower the side stability feet located toward the back on either side of the base plate until they touch the laboratory bench. Do not over-tighten to ensure level is maintained.
- 8 Reinstall the black caps on all of the screws.

Bath Temperature Probe Installation

- 1 Insert the water bath temperature probe into the BATH TEMP jack on the rear of the dissolution apparatus drive unit.



Figure 5. BATH TEMP Connector

- 2 Place the other end of the temperature probe through the hole in the center of the vessel plate in the water bath.

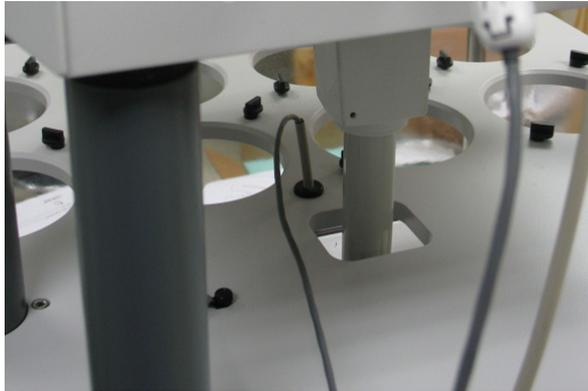


Figure 6. Positioning the Temperature Probe

Handheld Probe Installation

- 1 If applicable, insert the handheld vessel temperature probe cable into the VESSEL TEMP jack on the rear of the dissolution apparatus drive unit.
- 2 Place the vessel temperature probe in the hole in the vessel plate, at the left side rear of the apparatus.

Power Cord Connections

CAUTION

Ensure the correct voltage has been supplied prior to connecting the power cables. See “Environmental Requirements for Installation” on page 17.

- 1 Connect or verify connections of power cords to all components.
- 2 Plug the cords into electrical outlets of the appropriate voltage.

708-DS Initial Power Up

CAUTION

Ensure the retaining ties have been removed. Failure to do so could result in significant damage to the instrument.

- 1 Turn on the dissolution apparatus using the switch located on the left side of the drive unit.
- 2 Ensure the self-test conducted upon initialization passes, which is indicated by a green check mark.

Filling the Water Bath

- 1 If necessary, press and hold **Drive**  on the Main screen to allow for access to the water bath.
- 2 Fill the water bath to an appropriate level with purified water, using the label affixed to the bath as a guide. Be sure to allow for water displacement once the dissolution vessels are installed.

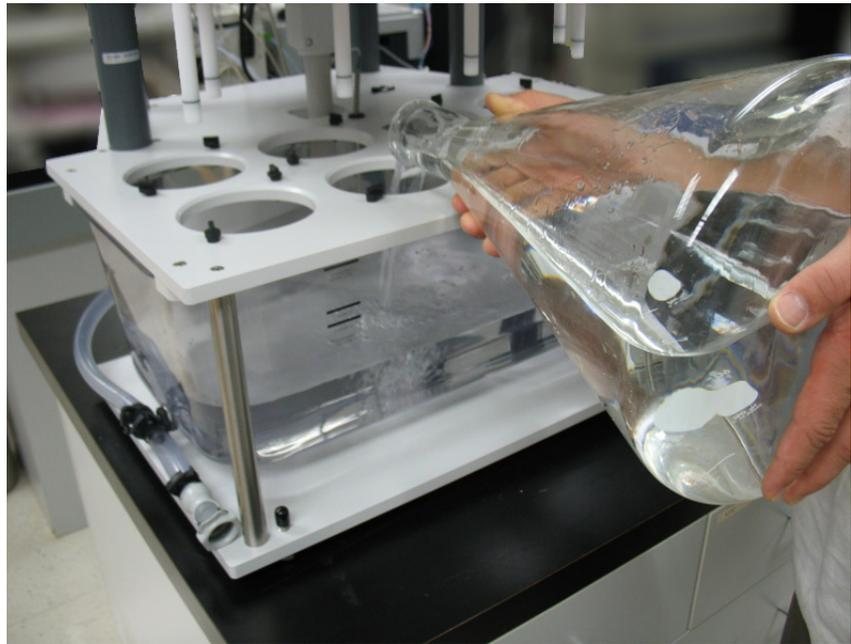


Figure 7. Filling the Water Bath

Level Instrument - Re-verification

- 1 With the water bath filled, re-verify the instrument level front to back and right to left using a bubble or digital level.

Heater / Circulator Power-up

NOTE

Use ultrapure water when possible to minimize scale and mineral buildup. Use algaecide to inhibit mold, bacteria, and algae growth. Check the label to ensure the formulation is compatible with the plastic materials used in the water bath construction. The flow paths of the heater / circulator are primarily stainless steel and should tolerate most clear water bath formulations.

- 1 Turn on the heater / circulator using the switch located on the unit.
- 2 Ensure water flow through the heater / circulator begins.
- 3 Ensure no leaks are present at any of the tubing connections.

Agilent 708-DS Setup

Complete the following sections to set up the Agilent 708-DS.

Vessel Installation

- 1 Press Drive  to move the drive unit to the fully raised position.
- 2 Carefully install the dissolution vessels in the vessel plate.
- 3 Engage the clips for each position over the rim of the vessel to secure it in place.
- 4 If necessary, place the blank vessel position covers on the vessel plate openings at Position 7 and Position 8. Rotate the two clips for each position over the cover to secure it in place.

Sample Manifold Installation

- 1 From the Main screen, press **Menu > Diagnostics**.

- 2 Position the sample manifold so the center screw is aligned with the center receptor beneath the drive unit. Ensure the left and right alignment shafts are also positioned correctly.

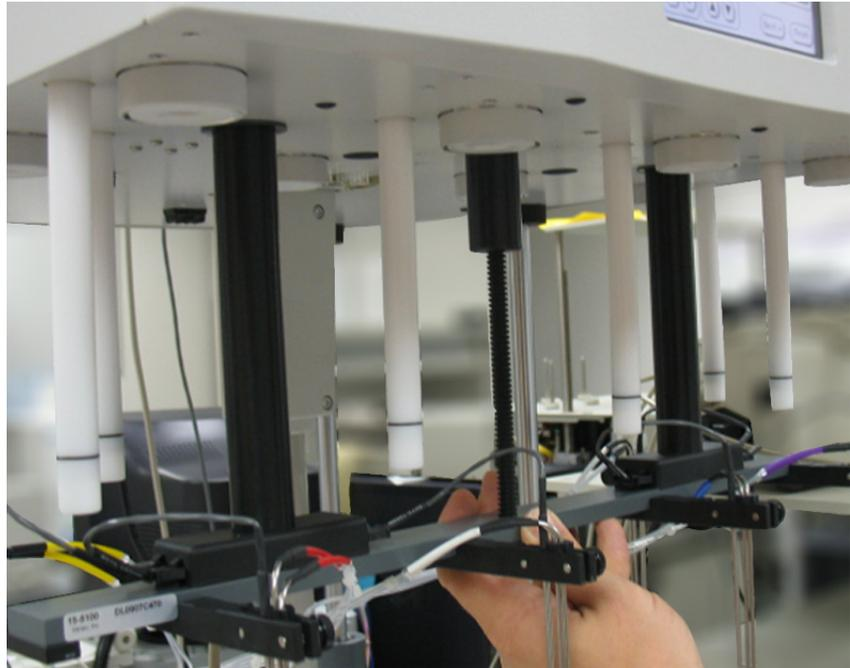


Figure 8. Positioning the Manifold for Installation

- 3 Push the manifold up into the drive unit until resistance is met.
- 4 While gently lifting upward, press Load from within the Manifold Control box. The internal drive should slowly lift the manifold to its fully raised position.

- 5 If the unit is equipped with AutoTemp, attach the two 9-pin RS232 cables to the underside of the drive unit.



Figure 9. Connecting AutoTemp

- 6 If the unit is equipped with autosampling, locate the black sample tubing fasteners and attach the fastener to the sample tubing wrap.

- 7 If applicable, affix the fasteners to the underside of the drive unit. A fastener is affixed to the left and to the right of the center drive unit lift.

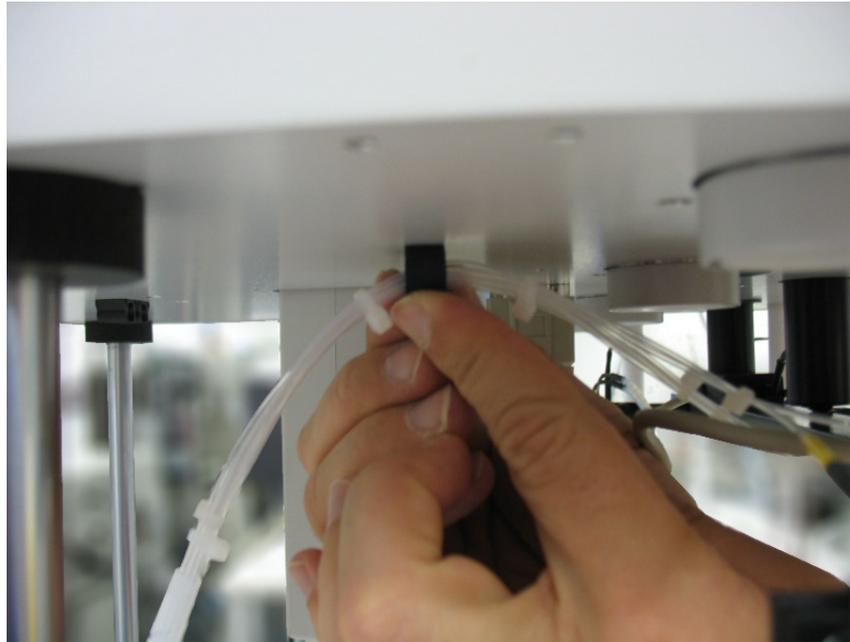


Figure 10. Affixing the Fasteners

Accessory Installation

- 1 Locate the appropriate accessories from the following list to be configured with the Agilent 708-DS: evaporation covers, Dosage Delivery Modules (DDMs), alignment posts, receptor shafts, paddle shafts, basket shafts, and shaft locking collars.
- 2 If necessary, press and hold **Drive**  to move the drive unit to its fully raised position.
- 3 If applicable, insert a receptor shaft (spin on/off shafts only) into each spindle location.
- 4 Slide an evaporation cover onto the shafts to be installed and insert the shaft into each spindle location.

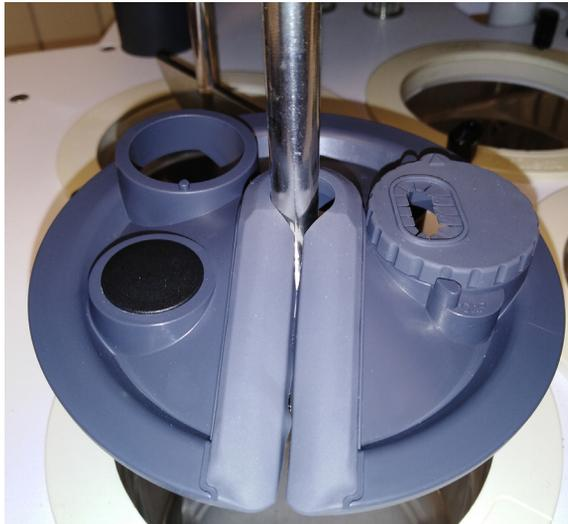


Figure 11. Installing Evaporation Covers on Shafts

If applicable, ensure the dial is rotated to the correct position for USP or ChP sampling. The cannulas and temperature probes should also be positioned accordingly on the sample manifold. This adjustment is made by shifting the individual arms of the manifold.

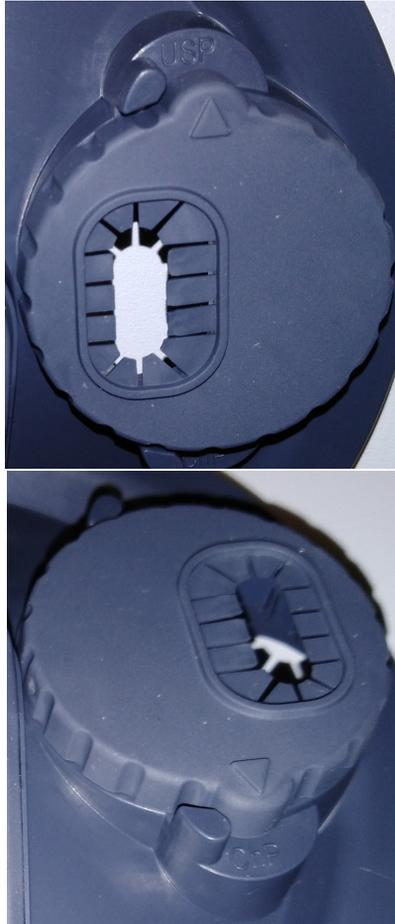


Figure 12. Evaporation Cover Dial (USP or ChP)

NOTE

Ensure that all sample tubing is behind the shaft and clear from all cables. See **Figure 13** below.

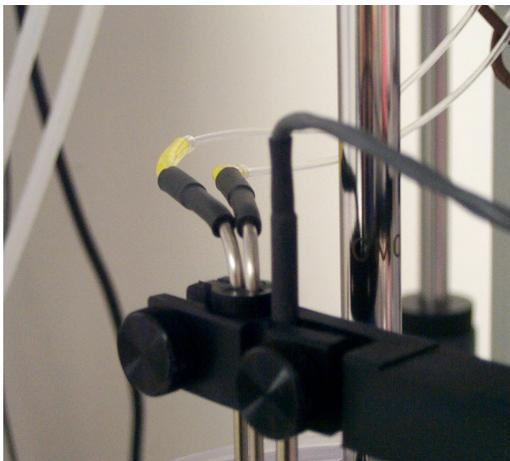


Figure 13. Proper Sample Tubing Position

- 5** Place a shaft locking collar on the top of each shaft and slide in down until it rests on top of the drive unit.



Figure 14. Shaft Locking Collars

- 6 Insert a DDM or alignment post into the DDM Alignment Shaft hanging from the underside of the drive unit.



Figure 15. Dosage Delivery Module and Alignment Post

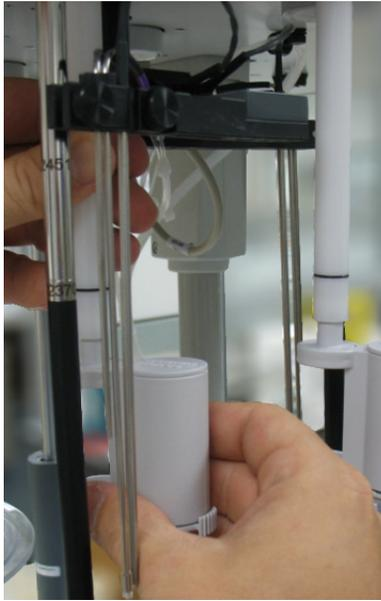


Figure 16. Inserting DDM into DDM Alignment Shaft

7 Repeat the previous step for all applicable positions.

- 8 Attach the DDM or alignment post to the evaporation cover by aligning the notch with the locking pin of the cover and pushing until fully seated.



Figure 17. Attaching DDM to Alignment Post

- 9 Repeat the previous step for all applicable positions.
- 10 Ensure that the sampling port of the evaporation cover is aligned properly with the sample manifold probes, if applicable. This may involve rotation of the dial to the appropriate USP or ChP sampling position and/or adjustment of the arms of the sampling manifold for 100/200 mL, 250 mL (CP), or 1-Liter vessels.

Setting Basket (Apparatus 1) Heights

- 1 Ensure the drive unit is fully raised and the basket shafts are pushed up sufficiently.
- 2 Clip the basket height gauge provided onto the bottom of the basket shaft.
- 3 Lower the drive unit to its operating position (until it stops).
- 4 With the shaft locking collar loosened, carefully lower the shaft until the bottom of the basket height gauge rests against the bottom of the vessel.
- 5 Ensure the shaft locking collar is flush against the top of the spindle assembly by rotating each shaft until resistance is met.
- 6 Tighten the shaft locking collar securely.



Figure 18. Tightening the Shaft Locking Collar

- 7 Raise the drive unit until sufficient clearance is available to move the basket height gauge to the next position.
- 8 Repeat steps 2 - 7 for all remaining positions.

Setting Paddle (Apparatus 2) Heights

- 1 Ensure the drive unit is fully raised and the paddle shafts are pushed up sufficiently.
- 2 Place a 25-mm height sphere in each vessel.

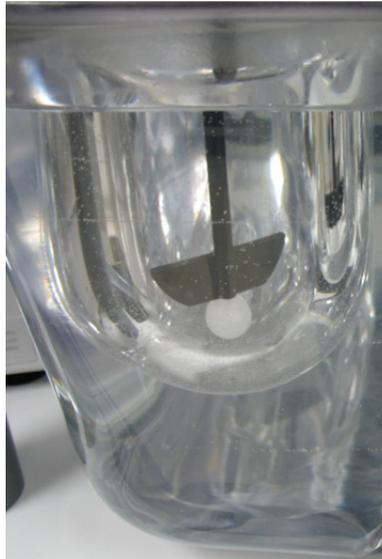


Figure 19. Inserting Height Spheres

- 3 Lower the drive unit to its operating position (until it stops).
- 4 With the shaft locking collars loosened, carefully lower each shaft until the bottom of the paddle blade rests against the height sphere.
- 5 Ensure the shaft locking collar is flush against the top of the spindle assembly by rotating each shaft until resistance is met.
- 6 Tighten each shaft locking collar securely.
- 7 Raise the drive unit to its home position.
- 8 Remove the height spheres from the vessels.



4

Operating the 708-DS

Main Screen - Options	40
Menu Screen - System Menu	58
General 708-DS Conventions	86
Agilent 708-DS Notifications	90

Main Screen - Options

The following sections detail the operations available from the 708-DS Main screen.

Main Screen - Display Parameters

The Main screen displays the system status and is updated once per second.

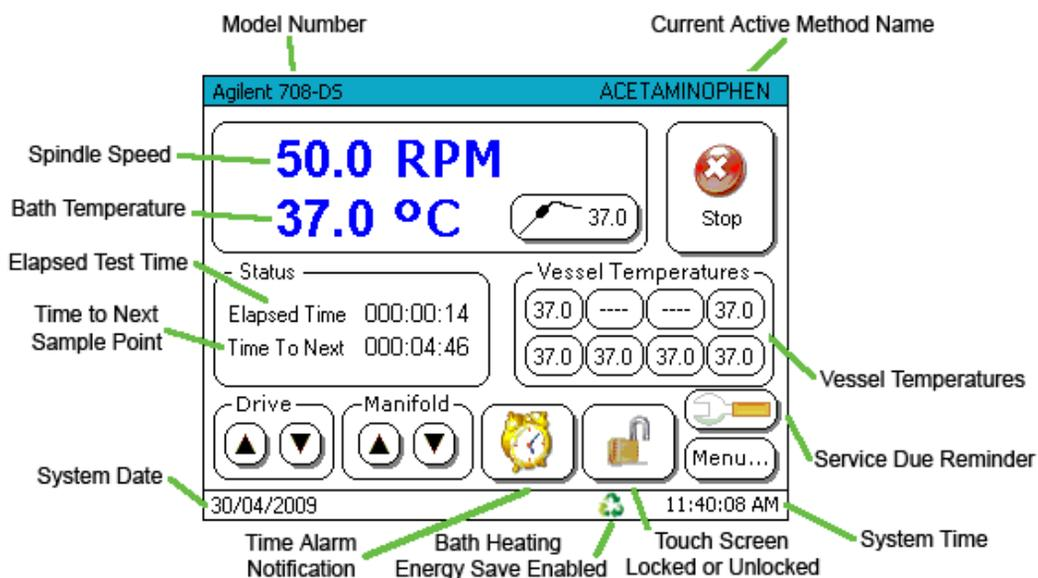


Figure 20. Main Screen - Display Parameters

NOTE

If the apparatus is in idle mode or running a manual method, press the area over the displayed RPM to quick-set the spindle speed. Upon pressing this area, the numeric data entry screen displays. Enter a valid number to change the RPM and press **Ok** to return to the Main screen. The spindle speed changes only if the spindle was previously running. If the spindle was idle, it starts.

Similarly, if you press the area over the displayed temperature, you can quick-set the bath temperature. Upon pressing this area, the numeric data entry screen displays. Enter a valid number to set the bath temperature setpoint. Press **Ok** to return to the Main screen.

Main Screen - Sample Temperatures

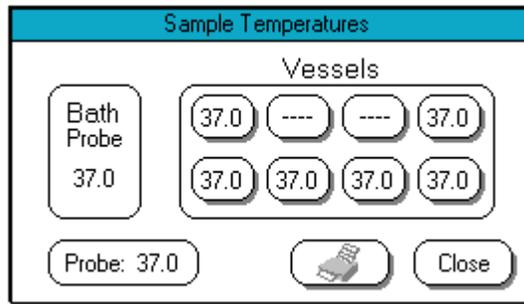


Figure 21. Sample Temperatures

NOTE

Sample Temperatures is available only if **Vessel Probe** is selected in the manufacturing configuration.

When you press , the Sample Temperatures box displays. The current bath temperature is displayed in the Bath Probe region and the current vessel probe temperature is displayed in the Probe region. Initially, the temperature at all vessel locations is displayed as "----".

Place the vessel probe in the appropriate vessel and press the corresponding vessel location on the screen to record the temperature of the vessel. If a vessel location is invalid (less than 8 active vessels), the unit beeps and the selected location remains blank.

NOTE

Repeatedly pressing a vessel location causes it to toggle between the current Probe temperature and "----". This allows you to correct or update the current temperature entry.

Press **Printer** to print the temperatures. If no printer is installed, the printer icon is not displayed. Press **Close** to close the window and return to the Main screen.

Main Screen - Run

When you press **Run**, the Select Start Type screen displays:

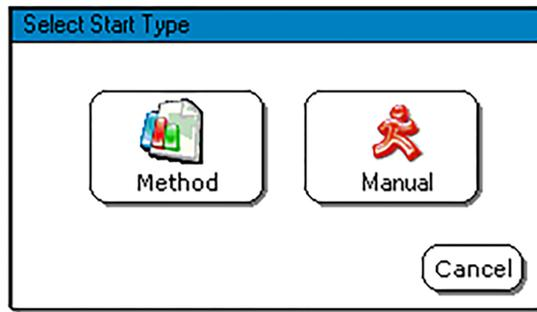
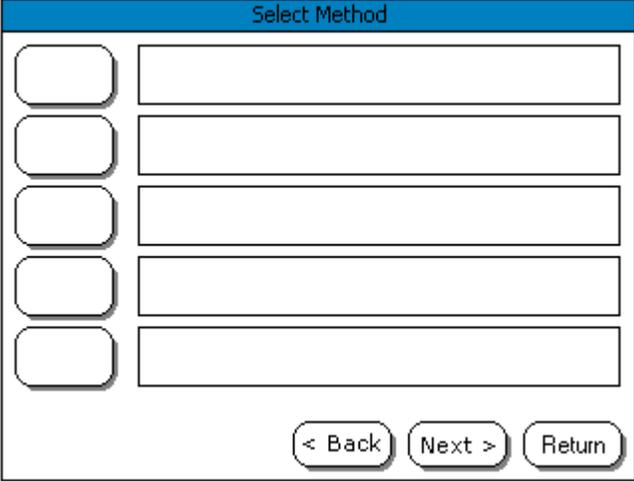


Figure 22. Select Start Type

From the Select Start Type screen:

- Press **Method** to select a method.
- Press **Manual** to start manual operation.
- Press **Cancel** to return to the Main screen.

Run - Select Method (Automated)

The screenshot shows a window titled "Select Method". Inside the window, there are five empty rectangular input fields arranged vertically. To the right of each input field is a small, empty rounded rectangular button. At the bottom right of the window, there are three larger rounded rectangular buttons labeled "< Back", "Next >", and "Return".

Figure 23. Select Method

You can enter up to 35 automated methods through the Select Method screen. In order to select a method to run, press the number of the method. Selecting a method takes you to the Start Options screen. Press **Return** to return to the Select Start Type screen.

Method - Start Options

Figure 24. Start Options

Option	Function
Instant	The method runs immediately after you press Ok .
Vessel Temperature	The manifold lowers the temperature probes into the media. When the selected vessel start temperature has been reached, the method starts. If no manifold is present, this option cannot be selected. For USP Apparatus 1 methods, you may install the 3-fin baskets to better equilibrate the media; the user will then be prompted to install the standard baskets with dosage forms to begin the test.
Bath Temperature	The method starts once the water bath has reached its selected start temperature.
Manual Dosage	Place a check in this box to indicate dosage forms will be introduced manually. A visual display will appear on-screen indicating when the user should drop dosage forms into each vessel.
Time Delayed	Displays the Date and Time screen. The apparatus begins the programmed test start sequence at the entered time and starts when the selected conditions have been reached.

From the Start Options screen:

- Press **Ok** to start the test with all operational parameters being monitored and return to the Main screen with the Stop button enabled.
- Press **Cancel** to abort the test start and return to the Select Method screen.

Run - Manual Operation

Figure 25. Manual Operation

Parameter	Range	Resolution
Temperature (Bath or Vessel)	25 - 55 °C	0.1 °C
Spindle Speed	10.0 - 250.0 RPM	0.1 RPM
Media Volume	100 to 2000 mL	Depends on configuration
(Test) Duration	000:00:10 to 999:59:59	1 second
Profile Print Interval	000:01:00 to 024:00:00	1 second
DDM Increment	000:00:00 to 001:39:59	1 second
Apparatus Type	1, 2, 5, or 6	
Option	Function	
Manual Dosage	Place a check in this box to indicate dosage forms will be introduced manually. A visual display will appear on screen indicating when you should drop dosage forms into each vessel.	

The Apparatus Type and Volume selection is used to control the manifold sampling depth when the manifold option is installed.

The Manual Operation screen allows you to run a manual method. Press the rectangular area following the text description to enter parameters. The Data Entry screen displays, allowing you to enter the data.

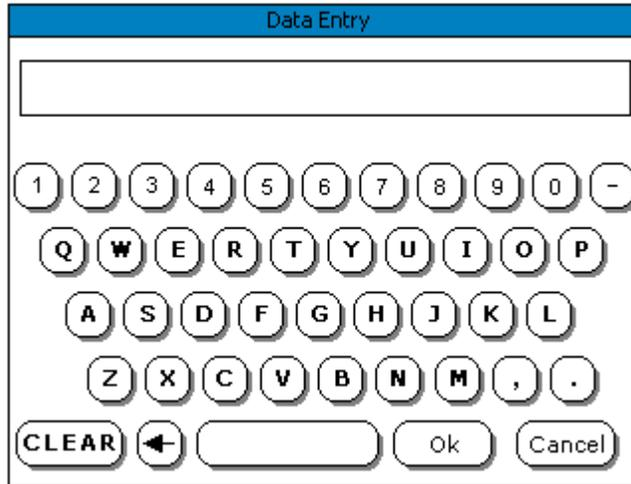


Figure 26. Data Entry

- Press **Cancel** to return to the Select Start Type screen without saving the test information.
- Press **Accept** to save the entered values and return to the Main screen.
- Press **Run** in the Manual Operation screen. The Main screen displays and the following actions take place:
 - a The Manually Insert Dosage(s) dialog box displays (“**Load Dosage Prompt**” on page 94).
 - b If DDMs are installed, tablets are automatically dropped into the vessels simultaneously or sequentially based on the DDM Increment. For example, a DDM Increment of 5 seconds drops the first tablet immediately, the second tablet 5 seconds after the first, and so on until all the tablets have been dropped. If the increment is zero, all tablets are dropped simultaneously.
 - c Elapsed time is reset to zero and incremented once per second while the test runs.
 - d The spindles start rotation at the set speed.

- e The bath temperature is maintained at the set temperature. If the bath is not at the set temperature when the test is started, the bath temperature is displayed in red while it heats or cools. A Bath Temperature Error ("**Bath Temperature Error**" on page 98) displays if the temperature is out of tolerance.
- f The displayed system date and time are updated once per second.
- g The measured rotational spindle speed and bath / vessel temperature(s) are displayed at an update rate of once per second.
- h Run is relabeled Stop.
- i If a printer is installed, test data is periodically printed based on the Profile Print Interval.
- j The test stops when the elapsed time equals the Test Duration.

Upon completion, a message is printed (if installed and enabled). The display reverts to the Main screen with the Run button enabled.

Main Screen - Stop

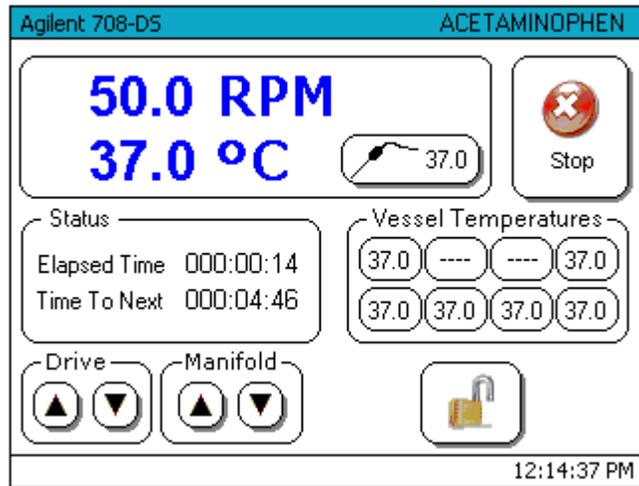


Figure 27. Main Screen - Stop

When the system is running, the Stop button is displayed. When you press **Stop**, the Pause Method screen displays:

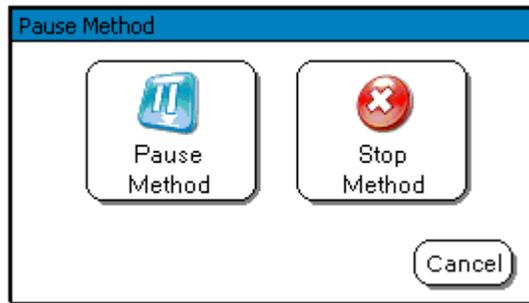


Figure 28. Pause Method

When this screen displays, the currently running method or manual test continues to run until you select an option. You have the option of pausing or stopping the test, or canceling this dialog and returning to the active method. The Pause Method button allows for media change during a dissolution test. If you select **Pause Method**, the Method Paused screen displays:

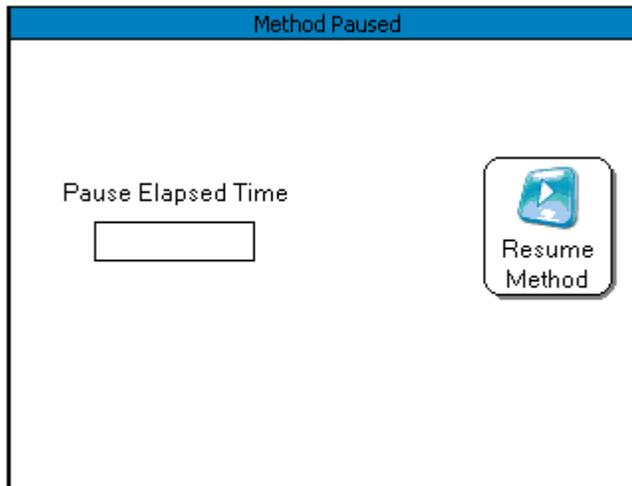


Figure 29. Method Paused

When a method is paused, the system takes the following actions:

- The spindles stop.
- If Full Media Change has been selected as a method parameter, the elapsed time counter stops counting.

- A Pause Elapsed Time counter starts counting. If the paused time exceeds the time specified in the Pause Duration Alarm, the Alarm warning screen displays. See “Pause Duration Exceeded” on page 98..

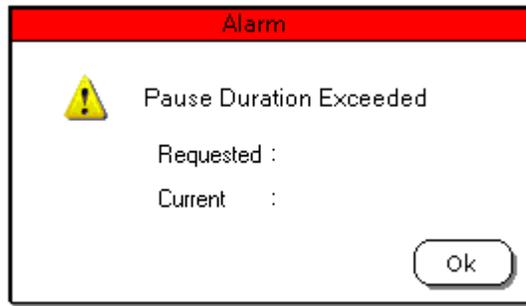


Figure 30. Pause Duration Alarm

A corresponding message is also printed for documentation. Clearing the Pause Alarm returns you to the Method Paused screen where you can complete the following tasks.

- The main drive may be raised or lowered.
- If installed, the manifold may be raised or lowered.
- If you press **Resume Method**, the system continues with the test. The paused time is printed, the Main screen displays, and the elapsed time counter resumes counting.

If you select **Stop Method**, the currently running test is terminated and the following actions are taken:

- The instrument returns to the Main screen.
- Stop is relabeled Run.
- The spindles stop.
- The elapsed time is reset to zero.
- The bath is maintained at the last entered temperature.
- The time, date, and the message TEST ABORTED BY USER are sent to the printer.

Main Screen - Drive Unit Up / Down

Press **Drive**  to raise the drive unit as long as the button is held. Release **Drive**  to stop the motion. Press **Drive**  to lower the drive unit. Double-tap **Drive**  to allow the drive unit to raise to its highest position without holding your finger on the button. There is no double-tap function for **Drive** .

NOTE

If the drive unit reaches the upper- or lower-limit switches, the motion stops automatically. Once an end-of-travel limit has been reached, only the arrow in the opposite direction operates.

For example, if the drive unit is raised completely, the Up arrow no longer causes motion but the Down arrow may be used to lower the drive unit. Once away from the limit sensors, both arrows resume normal operation.

Manual Sampling

If your Agilent 708-DS is not equipped with the Auto Sampling option for unattended sample collection, you will have to retrieve the samples manually. The evaporation cover design provides easy access to the appropriate dissolution sampling zone with the traditional manual sampling cannula.



Figure 31. Taking a Manual Sample

The 708-DS manual sampling bracket may also be used for manual sampling. This bracket is installed on the front of the unit and provides a repeatable, simplified mechanism for pulling samples. The bracket may be configured with the automated sampling manifold or used with resident probes.



Figure 32. Manual Sampling Bracket

Sampling and timing notifications are displayed at each sample timepoint based on the dosage introduction interval of the method. This display is accompanied by an audible alarm to alert the user to the upcoming event.

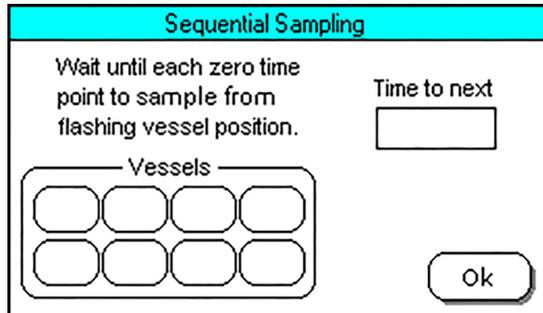


Figure 33. Sequential Sampling

Main Screen - Manifold Up / Down

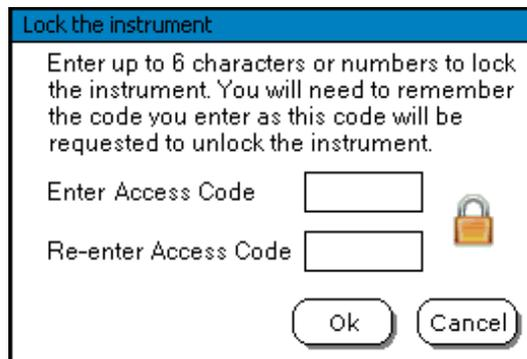
Press **Manifold**  to raise the manifold to its highest position. You do not need to hold the button to fully raise the manifold.

Press **Manifold**  to lower the manifold to the lowest position for the currently specified apparatus type. Setting the proper manifold depth for the appropriate volume and apparatus combinations is specified in **"Menu Screen - Calibration"** on page 76.

Main Screen - Lock

If Individual User Access mode is enabled, press  to lock the instrument.

If Individual User Access mode is disabled, press  to display the Lock the instrument screen.



Lock the instrument

Enter up to 6 characters or numbers to lock the instrument. You will need to remember the code you enter as this code will be requested to unlock the instrument.

Enter Access Code

Re-enter Access Code



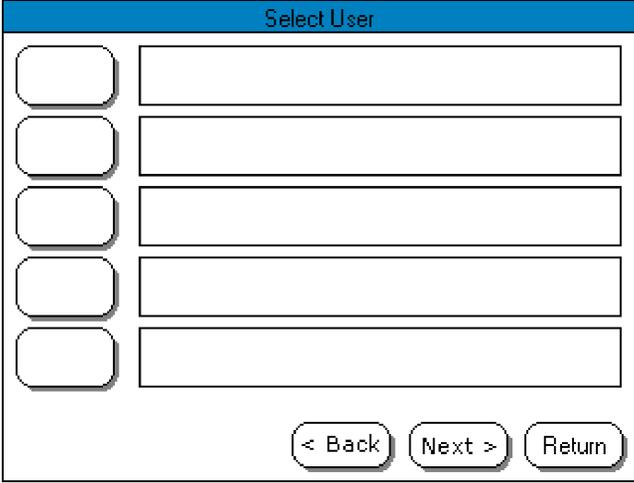
Figure 34. Lock the Instrument

Enter the access code, confirm it, and press **Ok**. The instrument returns to the Main screen and the Instrument Locked icon is displayed. Further changes to the instrument are prohibited until the instrument is unlocked or the power is cycled.

Main Screen - Unlock or Log In

Individual Access Mode Enabled

To unlock the instrument, press . The Select User screen displays.



The image shows a screen titled "Select User" with a blue header. Below the header, there are five rows, each consisting of a rounded rectangular button on the left and a rectangular text input field on the right. At the bottom right of the screen, there are three buttons: "< Back", "Next >", and "Return".

Figure 35. Select User Screen

Select the user to display the password screen. Enter the appropriate password and press **Ok**.

If the user is not yet created, an administrator must create the user. See "User Configuration" on page 84. for User Configuration creation details.

Individual Access Mode Disabled

The  icon indicates that the instrument is currently locked. In order to unlock the instrument, press . The Unlock the instrument screen displays requesting an access code:

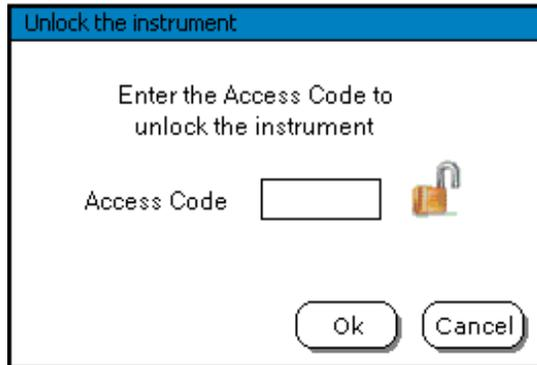


Figure 36. Unlock the Instrument

Enter the access code and press **Ok**. The instrument returns to the Main screen and  displays indicating that the system is unlocked. Full functionality of the instrument is now available through the front panel.

NOTE

If you forget your access code, the only way to regain access is to cycle the power.

Main Screen - Remote Control

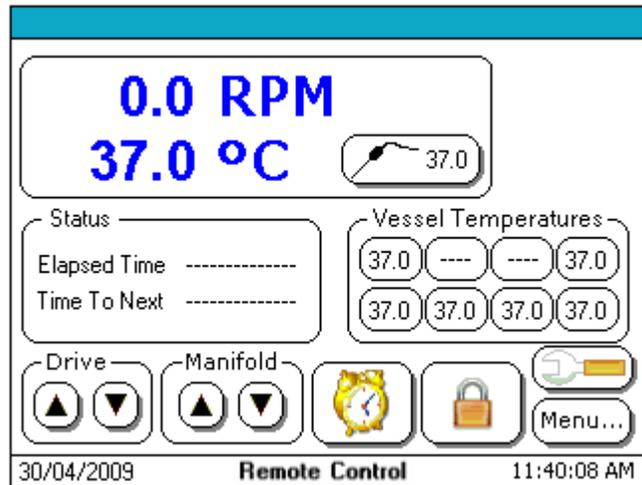


Figure 37. Remote Control

When the unit is being controlled remotely, the words Remote Control flash on the bottom of the screen, the Run / Stop button is not displayed, and the screen is locked. Although all other buttons are displayed, they are disabled.

Main Screen - Alarms

The alarm functions are accessed by pressing  on the Main screen. This notification applies only to the basic Timer Alarm function. The icon is displayed only if a Timer Alarm duration has been entered and the timer duration has not elapsed.

When the Timer Alarm duration elapses, an Alarm screen displays to notify the user and an alarm briefly sounds. The screen displays until you press **Ok**. See “Alarm Time Expired” on page 95..



Figure 38. Alarm Time Expired

NOTE

The Timer Alarm is provided as a convenience and does not affect the running of a manual test or an automated method. Other functions of the instrument may override the notification of the Timer Alarm if multiple events occur at the same time.

Main Screen - Maintenance Due Icon

When maintenance is due, a warning displays. Maintenance may include calibration or preventative Maintenance (“**Instrument Calibration Due**” on page 93). Once the warning has been acknowledged, the  icon displays as a reminder until the maintenance occurs. This icon is displayed on the Main screen until the appropriate due date has been updated under **Menu > Calibration**. See “Menu Screen - Calibration Screen 1” on page 76.

Menu Screen - System Menu



Figure 39. System Menu

Press **Menu** from the Main screen to access the System Menu. The System Menu provides an entry point for setting various system parameters. Each of the different functions available through this screen is described in detail in the following sections:

Method Editor	"Menu Screen - Method Editor" on page 59
Reports	"Menu Screen - Reports" on page 67
Alarms	"Menu Screen - Alarms" on page 70
Instrument	"Menu Screen - Instrument" on page 71
Calibration	"Menu Screen - Calibration" on page 76
Diagnostics	"Menu Screen - Diagnostics" on page 79
User Access	"Menu Screen - User Access" on page 81
Mfg. Config	"Menu Screen - Mfg. Config" on page 85

Menu Screen - Method Editor



The Method Editor allows you to create a new method, enter new method parameters, or modify an existing method.

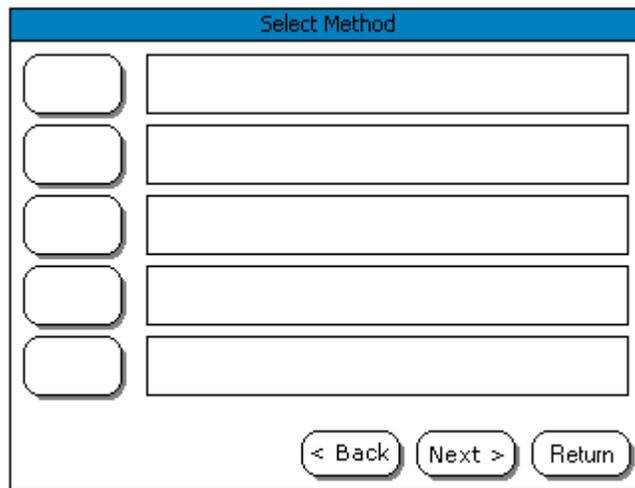


Figure 40. Method Editor

Thirty-five slots are allocated for storage of methods in system memory. In order to create a new method, select an empty memory slot, which is indicated by dashes in the name field. Tap an available method or slot. The Method Properties screen displays.

Method Properties Screen 1

Figure 41. Method Properties Screen 1

Option	Function
Name	32-character alphanumeric name used to describe the method.
Duration	The minimum time the method will take to run. The actual duration will be the value of this field or the sum of the final timepoint and final spin duration, whichever is greater.
Vessel Temp	Desired temperature of the vessels during a test.
Bath Temp	Desired temperature of the bath during a test.
Spindle Speed	Desired speed of the spindles during a test.
Media Volume	Volume of media in each vessel during a test.
Rotation Start Delay	Amount of time in seconds between the start of the test and the start of spindle rotation (maximum value of 10 seconds).
Apparatus Type	Currently installed apparatus.
Full Media Change	If this option is enabled, the elapsed time stops during a method pause. This allows for time to change the media. The elapsed time resumes counting when the method is resumed. If this option is not enabled, the elapsed time continues counting during a pause. If the Pause Duration Alarm time is exceeded, an alarm occurs.

From the Method Properties screen:

- Press **Clear > Yes** to return all method parameters to default values and return to the Select Method screen.
- Press **Cancel** to exit the routine without saving the values and return to the Select Method screen.
- Press **Save** to save the values to permanent memory and return to the Select Method screen.
- Press **Next** to save all values to permanent memory and proceed to the next Method Properties screen.

Method Properties Screen 2

Figure 42. Method Properties Screen 2

Option	Function
Enable Final Spin	Enables the final spin.
Final Spin RPM	Allows for spindle rotation at a speed different from the previous spindle speed.
Final Spin Duration	Duration of the final spindle speed setting.
DDM or Manual Increment	This is the time, in seconds, between successive tablet drops. Tablets are dropped starting with position one and continuing through all active vessels.
Profile Print Interval	This parameter controls the frequency that the spindle speed and temperature are printed. Entering a zero disables this function. Data is entered as HHH:MM:SS.

From the Method Properties screen:

- Press **Cancel** to exit the routine without saving the data and return to the Select Method screen.
- Press **Back** to save all values to permanent memory and return to the previous Method Properties screen.
- Press **Save** to save all values to permanent memory and return to the Select Method screen.
- Press **Next** to save all values to permanent memory and proceed to the next Method Properties screen.

Method Properties Screen 3

Figure 43. Method Properties Screen 3

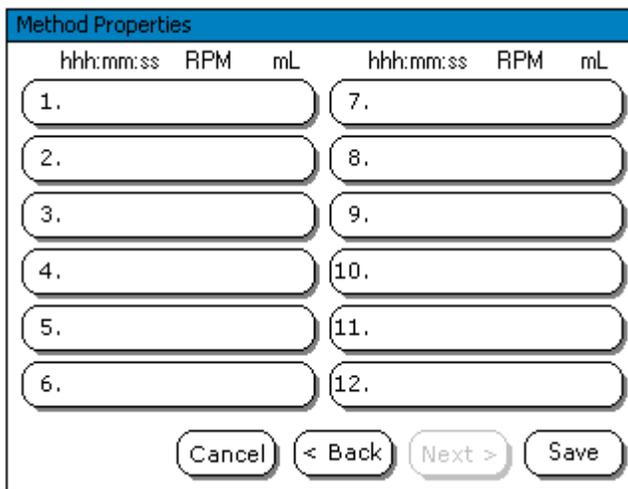
Option	Function
Enable Manifold	Enabling this option activates the optionally installed sampling manifold. The manifold lowers the sampling probes into the solution, allowing automated sample collection as well as temperature probes for automated temperature measurement.
Initial Temp	Enables recording and printing of the initial test temperature.
Final Temp	Enables recording and printing of the final test temperature.
Manifold Down Lead Time	This is the amount of time that the manifold lowers prior to the sample timepoint. This allows for equilibration of the temperature probes before vessel temperature measurement and for priming the system when an external pump is connected.

Option	Function
Manifold Down Duration	This is the amount of time that the manifold remains at its sampling position. The sum of the Lead Time and Duration must be greater than or equal to 30 seconds for accurate temperature readings.
Enable Sample Point Alarm	Enabling this option activates the sample notification dialog as well as activating the audible and visual alarms. This alarm draws the operator to the machine at the timepoint for manual sampling scenarios.
Alarm Lead Time	The Timepoint Approaching screen (" Timepoint Approaching " on page 91) is available to notify the user of upcoming timepoints. The time entered is how long before the timepoint the Timepoint Approaching screen displays. This alarm alerts the operator to an upcoming sample timepoint. Audible beeps will increase in frequency as the timepoint approaches.

From the Method Properties screen:

- Press **Cancel** to exit without saving the values and return to the Select Method screen.
- Press **Back** to save all values to permanent memory and return to the previous Method Properties screen.
- Press **Save** to save all values to permanent memory and return to the Select Method screen.
- Press **Next** to save all values to permanent memory and proceed to the next Method Properties screen.

Method Properties Screen 4

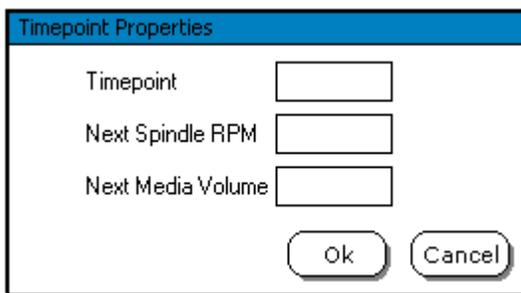


The Method Properties screen is a dialog box with a blue header. It contains two columns of input fields. The first column has six fields labeled 1. through 6., and the second column has six fields labeled 7. through 12.. Above the fields are two sets of column headers: 'hhh:mm:ss RPM mL' for the first column and 'hhh:mm:ss RPM mL' for the second column. At the bottom of the dialog are four buttons: 'Cancel', '< Back', 'Next >', and 'Save'.

Figure 44. Method Properties Screen 4

Using this Method Properties screen, you can specify up to twelve timepoints per test. These timepoints specify when changes in spindle speed, collection of samples, and recording of data occur.

Specify a timepoint by pressing the button for the corresponding timepoint. The Timepoint Properties screen displays requesting the new spindle speed and time for when that timepoint will be executed.



The Timepoint Properties dialog box has a blue header. It contains three input fields: 'Timepoint', 'Next Spindle RPM', and 'Next Media Volume'. At the bottom are two buttons: 'Ok' and 'Cancel'.

Figure 45. Timepoint Properties

You can enter a different RPM and vessel volume at each timepoint. The run report will print out the RPM that was running immediately before the timepoint. At the specified timepoint, the RPM will change (if applicable).

All timepoints are relative to the start of the test. A timepoint of 000:10:00 allows for a sample to be taken 10 minutes after the start of the test. If a second timepoint is entered as 000:25:00, the next sample point is programmed 25 minutes into the test.

From the Timepoint Properties screen:

- Press **Ok** to accept the timepoint information.
- Press **Cancel** to abort the process and return to the Method Properties - Timepoints screen.

NOTE

If a timepoint overlap is noted during timepoint entry, a warning icon displays in the lower left of the Method Properties - Timepoints screen. Timepoint overlap typically occurs when the manifold down lead time plus the manifold down duration time exceeds the timepoint interval. Press the warning icon to display the Timepoint Overlap Error screen ("**Timepoint Overlap Error**" on page 98). The warning icon remains on the screen as long as there are timepoint errors. A method can be saved even it contains timepoint overlap errors.

WARNING

Methods that contain overlap errors do not execute properly.

If sample point alarms have been enabled, timepoints are announced via the Sample Point Alarm screen.

From the Method Properties screen:

- Press **Cancel** to exit without saving the values and return to the Select Method screen.
- Press **Back** to save all values to permanent memory and return to the previous Method Properties screen.
- Press **Save** to save all values to permanent memory and return to the Select Method screen.

Method Cancel Dialog

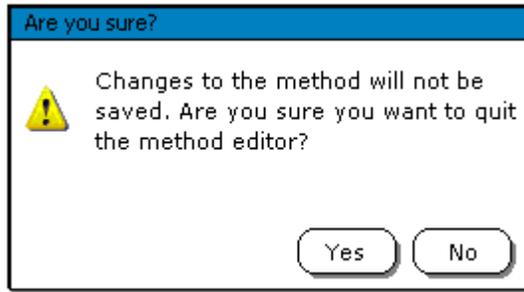


Figure 46. Are You Sure?

The Are you sure? screen displays when **Cancel** is pressed from one of the Method Properties screens. This option allows you to return to the Method Editor if Cancel is pressed inadvertently.

From the Are you sure? screen:

- Press **Yes** to return to the Select Method screen.
- Press **No** to return to the current Method Editor screen.

Menu Screen - Reports

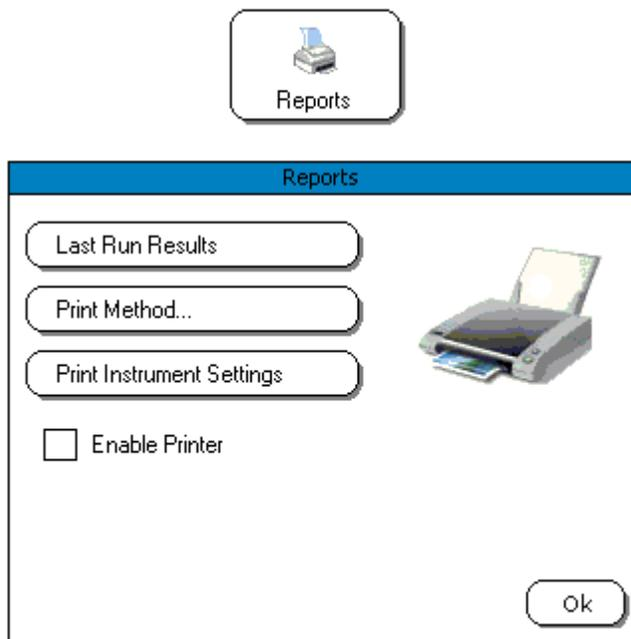


Figure 47. Reports

If a printer is installed, check **Enable Printer** to enable all printer functions. Uncheck the box to disable printing.

- To print results from the last completed test, press **Last Run Results**.

- To print the contents of a method, press **Print Method**. The Select Method screen displays prompting you to select a method to print:

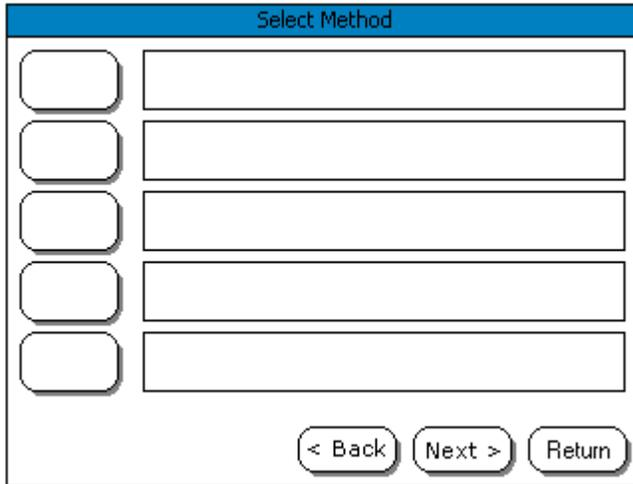


Figure 48. Select Method to Print

Available methods have a name associated with them. Methods that have not been configured have a series of dashes in the name field indicating available memory.

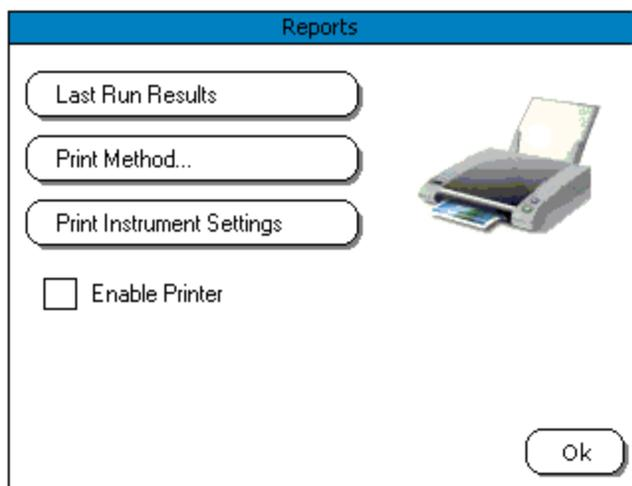


Figure 49. Reports

Press **Print Instrument Settings** to print all of the instrument configuration parameters and their current values. Press **Ok** to return to the Menu screen.

Menu Screen - Alarms

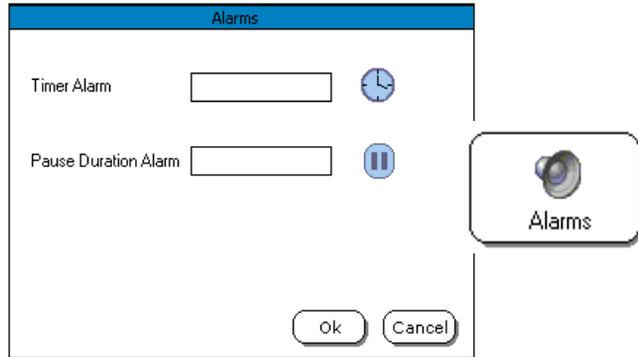


Figure 50. Alarms

Two different alarms are available through the Alarms screen:

Alarm	Function
Timer	This alarm functions as a countdown timer. Enter the duration as HH:MM:SS. Press Ok to start the timer. When the alarm duration expires, the Alarm Time Expired warning displays (" Alarm Time Expired " on page 95) and an audible alarm sounds. This alarm sounds until you press Ok . To cancel a previously running alarm before time has expired, enter 0 for the Timer Alarm value.
Pause Duration	This field allows you to set the maximum pause period while running a method. When a running method is paused, a timer starts. If the pause time exceeds this Pause Duration, the Pause Duration Exceeded error displays (" Pause Duration Exceeded " on page 98), an alarm sounds, and a message is sent to the printer. Entering a zero in this field disables the Pause Duration Alarm.

From the Alarms screen:

- Press **Ok** to enable the Timer Alarm and return to the Menu screen.
- Press **Cancel** to discard any changes and return to the Menu screen.

Menu Screen - Instrument



Menu Screen - Instrument Settings Screen 1

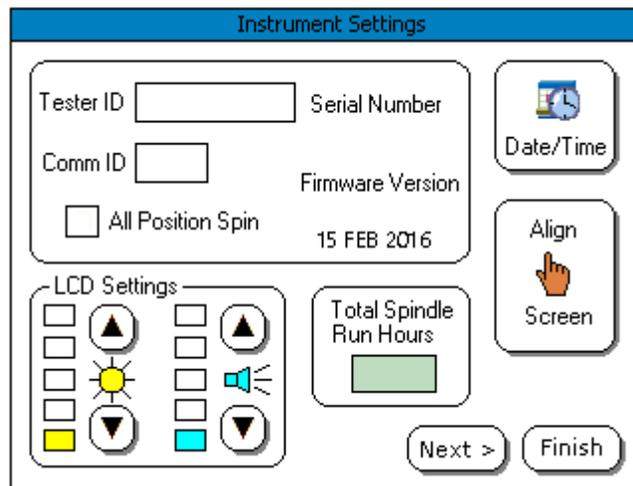


Figure 51. Menu Screen - Instrument Settings Screen 1

Option	Function
Tester ID	A twelve-character field that helps identify the instrument. The Tester ID is printed out with the other system information during a test.
Comm ID	The address of the instrument on the communications bus. This address is used by external host programs to communicate with the instrument. The Comm ID may range from 01 to 99.
All Position Spin	Allows the spindle to rotate regardless of the position of the drive unit. This is useful when calibrating spindle speed as the spindle may rotate while the drive unit is in its upper position. If this option is not checked, the spindle only rotates when the drive unit is in its lowest position.

Option	Function
Date / Time	Allows you to set the current date and time as well as set the desired format. Press the date format button to advance the date format to the next available format. Date formats are US (MM/DD/YYYY), ISO (YYYY/MM/DD) or Metric (DD/MM/YYYY). Time formats are either 12-hour (2:42:00 PM) or 24-hour (14:42:00).
Align Screen	Starts the touchscreen calibration routine. The screen prompts you to press three different circles in succession to calculate the calibration. To verify calibration, press the 4 circles in the screen corners. Multiple Xs normally display in each circle. Touchscreen calibration values are stored in permanent memory for further use. If the touchscreen calibration was entered inadvertently, do not touch the screen. After a 5-second delay, the calibration routine aborts and the touchscreen calibration is not changed.
LCD Settings	Sets the screen brightness and the audio feedback volume. Pressing either Up arrow increases the desired value. The corresponding bar graph increases to reflect the new setting. Pressing the Down arrow decreases the values.

From the Instrument Settings screen:

- Press **Finish** to return to the Menu screen. All values are stored in permanent memory.
- Press **Next** to advance to the next Instrument Settings screen.

Menu Screen - Instrument Settings Screen 2

The screenshot shows the 'Instrument Settings' screen. It has a blue header with the text 'Instrument Settings'. Below the header, there are two main sections: 'Tolerances' and 'Vessels'.
 - The 'Tolerances' section contains two rows: 'Temperature Tolerance' with an input box and '°C' to its right, and 'Speed Tolerance' with an input box and 'RPM' to its right.
 - The 'Vessels' section is split into two columns. The left column has radio buttons for '2000ml', '1000ml', '200ml', '100/250ml', '100ml', and '250ml'. To the right of these is the text 'Number Active' followed by an input box. The right column has two checkboxes: 'Enable Alarms and Errors' and 'Enable DDM'.
 - At the bottom of the screen, there are three buttons: '< Back', 'Next >', and 'Finish'.

Figure 52. Menu Screen - Instrument Settings Screen 2

Option	Function
Temperature Tolerances	Allowable temperature variation from the set point. When alarms are enabled in Instrument Settings, deviations greater than this will cause an alarm but not stop the method.
Speed Tolerance	Allowable spindle speed variation from the set point. When alarms are enabled in Instrument Settings, deviations greater than this cause an alarm but do not stop the method.
Vessels	This parameter sets the current installed vessel size (2000, 1000, 250, 200, or 100 mL). This value is used to determine various settings throughout the operation of the instrument but most importantly sampling manifold depth which can be programmed from the Calibration screen.
Number Active	Sets the number of vessels used in a test. The remainder of the vessels are ignored until this parameter is changed.
Enable Alarms	If checked, all alarms are activated. If not checked, all alarms are suppressed on the user interface and on a printout if a printer is installed.
Enable DDM	Enables the Dosage Delivery Modules. This setting overrides the method setting. If DDM is not enabled, the DDM Increment box is grayed out in Method Editor.

From the Instrument Settings screen:

- Press **Back** to return to the previous Instrument Settings screen.
- Press **Finish** to save all settings and return to the Menu screen.
- Press **Next** to advance to the next Instrument Settings screen.

Menu Screen - Instrument Settings Screen 3

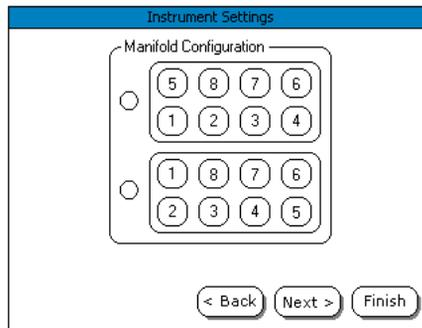


Figure 53. Menu Screen - Instrument Settings Screen 3

Option	Function
Manifold Configuration	Selects the numbering scheme to be used throughout testing.

From the Instrument Settings screen:

- Press **Back** to return to the previous Instrument Settings screen.
- Press **Finish** to save all settings and return to the Menu screen.

Menu Screen - Instrument Settings Screen 4

Figure 54. Menu Screen - Instrument Settings Screen 4

Option	Function
Bath Heating Energy Save	Enable this option to program the power-save feature for the heater/circulator. Heater/circulator on and off times can be specified to conserve energy and reduce stress on the components of the unit.

From the Instrument Settings screen:

- Press **Back** to return to the previous Instrument Settings screen.
- Press **Finish** to save all settings and return to the Menu screen.

Menu Screen - Calibration



Menu Screen - Calibration Screen 1

Paddles/Baskets Sampling Depth

Paddles

Baskets

Current Position

Manifold

Figure 55. Paddles/Baskets Sampling Depth

The Calibration screen allows you to fine-tune the height of the manifold for Apparatus 1 (Baskets) and Apparatus 2 (Paddles). The volumes shown change based on the installed vessel size set in Instrument Setting Screen 2 (“**Menu Screen - Instrument Settings Screen 2**” on page 72). The following vessel size ranges are configurable:

Size	Max	Min
2L	2000 mL	1650 mL
1L	900 mL	500 mL
250 mL	300 mL	100 mL
200 mL	200 mL	150 mL
100 mL	100 mL	80 mL

Each of the gray boxes contains a numeric value. The value is the sampling location of the manifold as measured from its home position. These values range from 0 at home to 999 at the manifold's lowest position. If no values have ever been entered, the default values, protected in the firmware, are shown for each Apparatus. If previous values have been entered, these are displayed. Default values for a 1L system are shown in the screen shot.

NOTE

When changing between the 100ml and 250ml options an “Are you sure?” dialogue is displayed as the saved calibration depths and timepoint volumes will be reset to default for the selected volume.

NOTE

When changing between the 100 mL and 250 mL options, the previously stored calibration depths and timepoint volumes will be reset to default for the selected volume.

To set a new position, press **Manifold** ▲ or **Manifold** ▼ to raise or lower the manifold. The manifold will move as long as the button is held. The current location of the manifold is displayed above **Manifold** ▲▼. Once the desired position is reached, press Save on the corresponding field. This updates the display for that field.

From the Paddles/Baskets Sampling Depth screen:

- Press **Restore Defaults** to replace all of the displayed values with the default values, which are protected in the firmware.

- Press **Cancel** to exit the routine without saving the values and return to the Menu screen.
- Press **Finish** to save all values to permanent memory and return to the Menu screen.
- Press **Next** to save all values to permanent memory and proceed to the next Calibration screen.

Menu Screen - Calibration Screen 2

The image shows a screen titled "Regulatory Calibration Date". It contains two text input fields: "Calibration Due Date" and "PM Due Date". To the right of these fields is a small calendar icon. At the bottom of the screen, there are four buttons: "Cancel", "< Back", "Next >", and "Finish".

Figure 56. Regulatory Calibration Date

Calibration and Preventative Maintenance due dates may be entered on the Regulatory Calibration Date screen. A reminder displays over the Main screen to indicate that calibration or maintenance is required. You must develop your own Calibration and Preventative Maintenance schedules.

From the Regulatory Calibration Date screen:

- Press **Cancel** to exit the routine without saving the values and return to the Menu screen.
- Press **Back** to save all values to permanent memory and return to the previous Calibration screen.
- Press **Finish** to save all values to permanent memory and return to the Menu screen.

Menu Screen - Diagnostics



Menu Screen - Diagnostics Screen 1

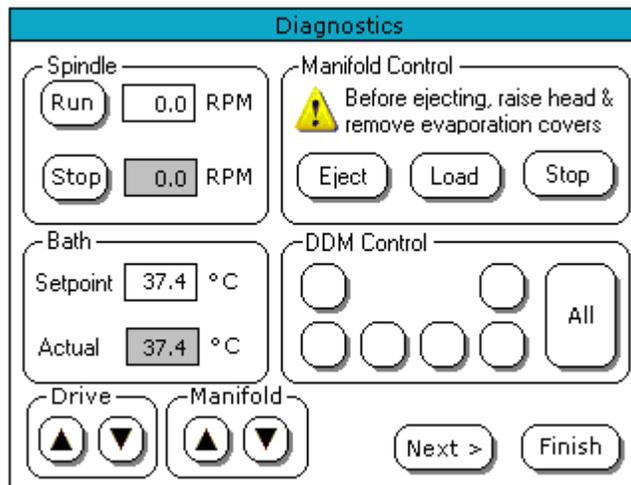


Figure 57. Diagnostics Screen 1

Option	Function
Spindle	Allows control of the spindle. Press the box next to Run and enter the desired RPM. Press Run to start the spindle and press Stop to stop the spindle. The actual RPM displays in the gray box below the RPM set point.
Bath	Allows control over the bath temperature. Press the box next to Setpoint and enter the desired temperature in the open field that displays. Press Ok to activate the temperature controller until the set point temperature is reached. The actual temperature is displayed in the gray box below the set point temperature.
Manifold Control	The manifold may be ejected for cleaning and maintenance. Before ejecting, raise the head, disconnect the two thermistor cables, remove the tubing from the tubing clamps, and remove the evaporation covers. Press Eject to lower the manifold until it is free. To reattach the manifold, insert it into the guides and gently lift it up while pressing Load. Release the manifold once it begins to rise on its own. Press Stop to stop all manifold motion.

Option	Function
Dosage Delivery Module Control	The DDMs are solenoid activated doors that hold a tablet above each vessel. Press the corresponding button to open a DDM. Press All to open all DDMs simultaneously.
Drive	Press Drive ▲ to raise the drive unit. Keep pressing it to raise the drive unit to the maximum height. Release the button to stop the motion. Press Drive ▼ to lower the drive unit.
Manifold	Press Manifold ▲ to raise the manifold to the maximum height. Press Manifold ▼ to lower the manifold to the appropriate position based on the apparatus type and volume.

From the Diagnostics screen:

- Press **Next** to proceed to the second Diagnostics screen.
- Press **Finish** to return to the Menu screen.

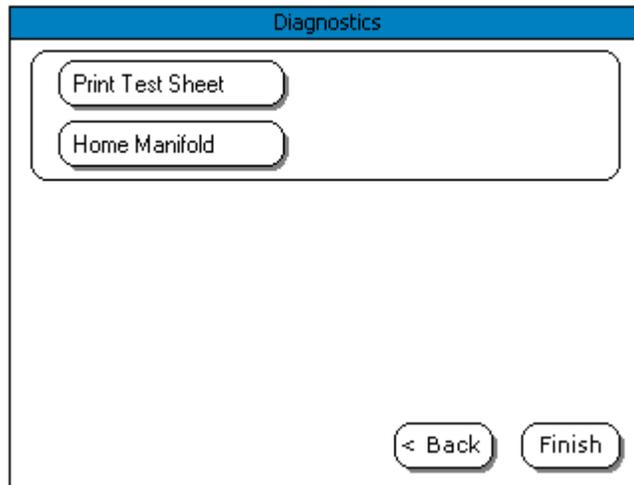


Figure 58. Menu Screen - Diagnostics Screen 2

Option	Function
Print Test Sheet	Sends a test message to the printer. A printer is required for this option.
Home Manifold	Automatically returns the sample manifold to its home (fully raised) position.

From the Diagnostics screen:

- Press **Back** to return to the previous Diagnostics screen.
- Press **Finish** to return to the Menu screen.

Menu Screen - User Access

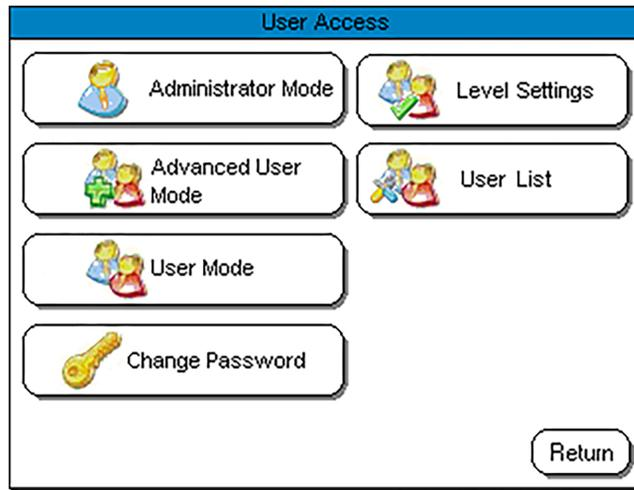


Figure 59. User Access

The User Access screen provides an Administrator with the ability to change the security settings of the 708-DS.

Option	Function
Administrator Mode	Provides access to system security settings and all functions enabled in column 1 of the User Level Settings (“ User Level Settings ” on page 83).
Advanced User Mode	Provides access to functions enabled in column 2 of the User Level Settings (“ User Level Settings ” on page 83).

Option	Function
User Mode	Provides access to functions enabled in column 3 of the User Level Settings (“ User Level Settings ” on page 83). Note: Instrument Mode Settings can be modified only by a user with the appropriate User Level. Only an Administrator can access certain security settings.
Change Password	Allows an Administrator to change the 708-DS password.

Option	Function
--------	----------

Level Settings

User Level Settings			
Function	1	2	3
Methods: Create / Edit / Delete	✓		
Lock Screen	✓	✓	✓
Alarms: Timer & Pause Duration	✓	✓	✓
Due Dates: Calibration & PM	✓	✓	
Diagnostics	✓	✓	
Instrument: LCD, Date/Time, Align Screen	✓	✓	✓
Instrument: Comm ID, All Pos. Spin, Man. Config.	✓	✓	
Instrument: Tolerances, Vessels, Enables	✓	✓	
Instrument: Energy Save	✓	✓	

1. Administrator, 2. Advanced User, 3. User

Ok

Figure 60. User Level Settings

You can use user levels to restrict changes to specific instrument settings and method parameters. Three levels of access are provided:

- Administrator
- Advanced User
- User

Option	Function
User List	Allows an Administrator to configure specific users for the 708-DS. Once a position is selected from the User List screen, the User Configuration screen displays.

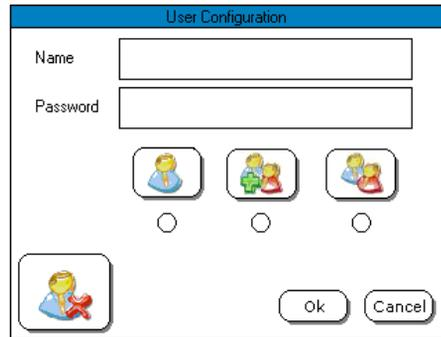


Figure 61. User Configuration

The User Configuration screen can be used to create or edit users. The screen contains settings for Name, Password, and User Level as well as the ability to delete the user.

If the Username/Password feature is enabled, operation of the 708-DS will be disabled from the Main screen until a valid user (or Administrator) logs in from the Main screen. A locked instrument can be identified by the  icon displayed on the Main Screen. The User Login screen displays when  is pressed.

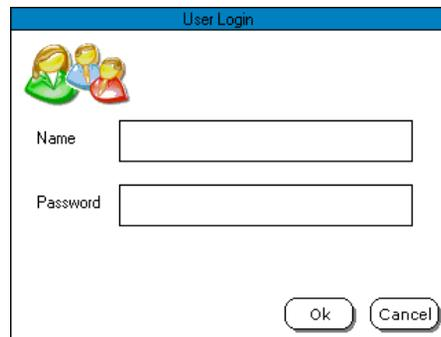


Figure 62. User Login

NOTE

You can customize each user access level, but changes can only be made when the instrument is in Administrator mode. The function settings of each user level are applied to instrument operation with and without the Username/Password feature enabled.

From the User Level Settings screen:

- Press **Ok** to return to the previous window.

Menu Screen - Mfg. Config

Password access is required to access the Manufacturing Configuration settings. Instrument options configured here include:

- Dosage delivery module
- Auto sampling
- Auto temp
- Username / Password Access
- Printer
- Vessel probe
- Language option
- Maximum number of vessels
- Administrator password reset

Contact Agilent Customer Care (“[Obtaining Warranty and Other Services](#)” on page 116) if access to these options is required.

General 708-DS Conventions

Alphanumeric Data Entry

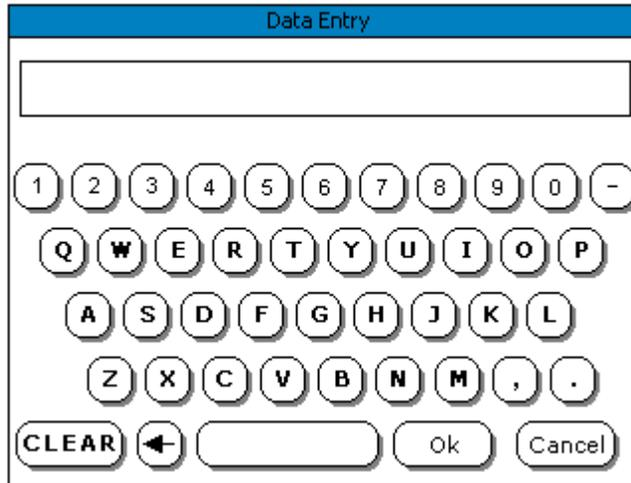


Figure 63. Data Entry

The Data Entry screen acts as on-screen keyboard for the instrument. The left arrow functions as a backspace key. Spaces, periods, and dashes are provided as usable characters.

From the Data Entry screen:

- Press **CLEAR** to remove all characters from the open field.
- Press **Ok** to return to the previous window. The data entered is returned and placed in the appropriate data entry box.
- Press **Cancel** to clear the entered data and return to the previous window.

Numeric and Time / Date Entry

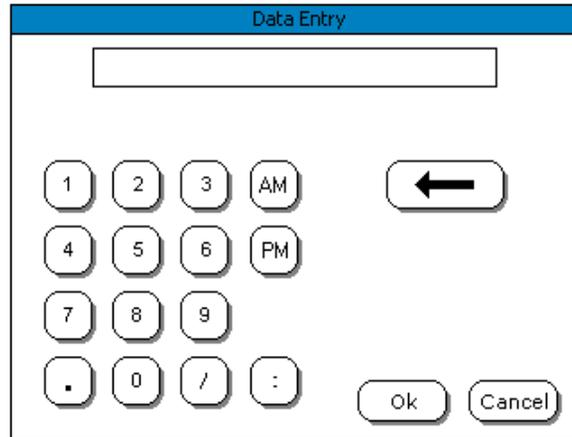


Figure 64. Time / Date Entry

This screen is a specialized subset of the alphanumeric data entry screen. It has been optimized for numeric, date, and time data entry. As keys are pressed, they display in the upper box of the display. The valid data entry range and the previous data entry will be shown just below the display. The maximum valid data string length is 25 characters. Special function keys are defined below:

Option	Function
←	Backspaces and deletes one character every time it is pressed.
/	Used as separator in date entries.
:	Used as separator in time entries.
AM	Designates morning when entering time in 12-hr format.
PM	Designates afternoon when entering time in 12-hr format.

From the Numeric and Time / Date Entry screen:

- Press **Ok** to return to the previous window. The data entered is returned and placed in the appropriate data entry box.

Parameter Limits / Ranges

Parameter	Range	Resolution	Configured
Spindle Speed	10.0 - 250.0 RPM	0.1 RPM	"Main Screen - Display Parameters" on page 40 "Run - Manual Operation" on page 45 "Method Properties Screen 1" on page 60 "Menu Screen - Diagnostics Screen 1" on page 79
Temperature (Bath or Vessel)	5 - 55 °C	0.1 °C	"Main Screen - Display Parameters" on page 40 "Method Properties Screen 1" on page 60 "Menu Screen - Diagnostics Screen 1" on page 79
Serial Number	15 characters		Factory
Profile Print Interval	000:01:00 to 024:00:00	1 second	"Run - Manual Operation" on page 45 "Method Properties Screen 2" on page 61
Test Duration	000:00:10 to 999:59:59	1 second	"Run - Manual Operation" on page 45 "Method Properties Screen 3" on page 62
Final Spin Duration	000:00:10 to 001:00:00	1 second	"Method Properties Screen 2" on page 61
Tester ID	12 characters		"Menu Screen - Instrument Settings Screen 1" on page 71
Comm ID	01 to 99	01	"Menu Screen - Instrument Settings Screen 1" on page 71
Media Volume	100 to 2000 mL	1 mL	"Run - Manual Operation" on page 45 "Method Properties Screen 1" on page 60

Parameter	Range	Resolution	Configured
DDM Increment	000:00:00 to 001:39:59	1 second	"Run - Manual Operation" on page 45 "Method Properties Screen 2" on page 61
Temp Error Tolerance	0 to ± 0.5 °C	1 second	"Menu Screen - Instrument Settings Screen 2" on page 72 Note: A value of 0 disables the tolerance check.
Speed Error Tolerance	0 to 1.0 RPM	1 second	"Menu Screen - Instrument Settings Screen 2" on page 72 Note: A value of 0 disables the tolerance check.
Manifold Down Time	000:00:00 to 001:00:00	1 second	"Method Properties Screen 3" on page 62
Manifold Lead Time	000:00:00 to 000:01:00	1 second	"Method Properties Screen 3" on page 62
Sample Point Alarm	000:00:05 to 000:02:30	1 second	"Method Properties Screen 3" on page 62
Method Pause Time	00:00 to 1:00:00	1 second	"Menu Screen - Alarms" on page 70
Timer Alarm	00:00 to 24:00:00	1 second	"Menu Screen - Alarms" on page 70

Agilent 708-DS Notifications

Main Screen - Error Conditions

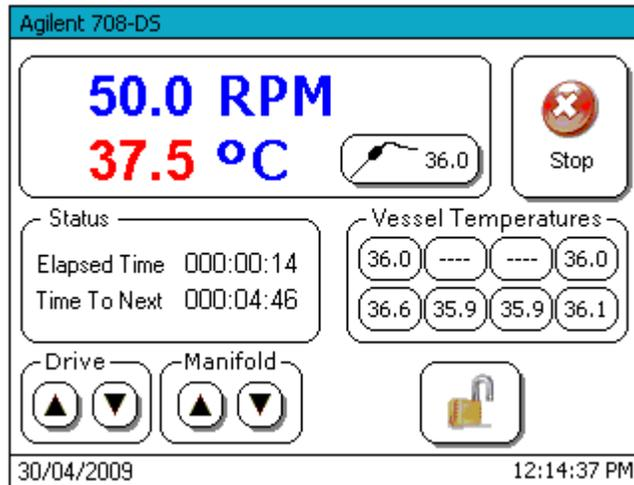


Figure 65. Main Screen with Red Error Conditions

Tolerance limit monitoring is enabled when the spindle starts moving and the bath starts heating. Any fluctuations beyond the tolerance limits cause the corresponding displayed value to turn from blue to red and a dialog box with the corresponding error message to be displayed.

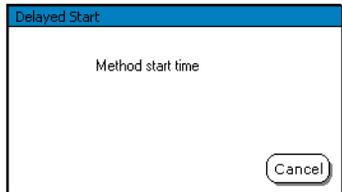
More critical errors are displayed on a separate error screen. These errors may include the following:

- Bath temperature out of tolerance
- Bath probe open circuit
- Bath probe short circuit
- Heater probe open circuit
- Heater probe short circuit
- Heater to bath probe offset > 6°. For this error, it is important to disable the heater to prevent potential damage due to a dry system.

- Spindle speed error if the set speed and actual speed differ by a value greater than the spindle speed tolerance.

If an error occurs during a test, you must clear the error screen manually. Clearing the error screen also clears the audible alarm. If a printer is installed and alarms are enabled in the Instrument Settings, the error condition is printed for record keeping.

Dialog Screens (Green / Blue)



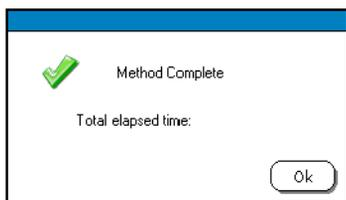
Delayed Start Dialog

The Delayed Start screen displays when a delayed start time has been entered (“**Method - Start Options**” on page 44). The date and time when the method starts is displayed on the screen. Access to the remainder of the instrument functions is prohibited until the method has finished. Press **Cancel** to stop the loaded method and return to the Main screen.



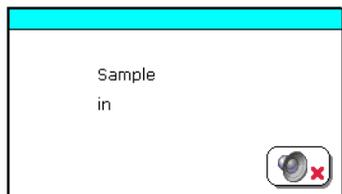
Delayed Start Date / Time Dialog

Displays when Time Delayed Start has been selected (“**Method - Start Options**” on page 44). Press the box beside either the Date or Time prompts to display the Numeric and Date and Time Entry screen (“**Numeric and Time / Date Entry**” on page 87). Press **Cancel** to abort the process and return to the previous screen.



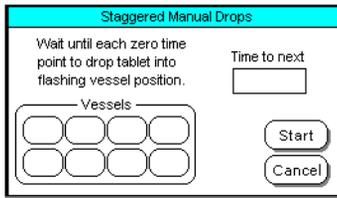
Method Completion

The Method Completion screen displays when a method has completed. The total elapsed time, current date, and time the method completed is displayed. Press **Ok** to return to the Main screen.



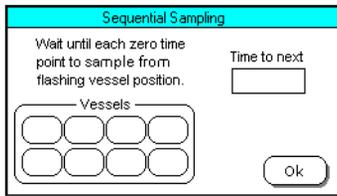
Timepoint Approaching

Displays when a sample point is approaching (“**Method Properties Screen 4**” on page 64). The time to the next sample point is indicated on the screen. Press **Ok** to clear the dialog and return to the Main screen. If no action is taken, this screen clears itself when 0 seconds is reached.



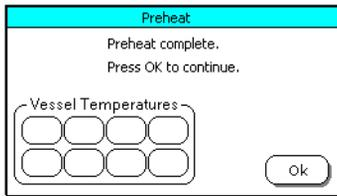
Staggered Manual Drops

The Staggered Manual Drops screen provides a visual indication of when to drop tablets for manual introduction.



Sequential Sampling

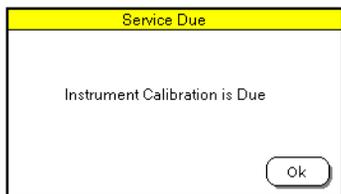
The Sequential Sampling screen provides visual indication of when to remove sample manually from each dissolution vessel. This interval is automatically determined based on the dosage introduction interval of the method.



Preheat Complete

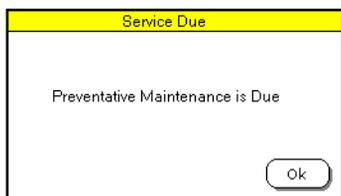
Displays when the preheat process has completed.

Warning Screens (Yellow)



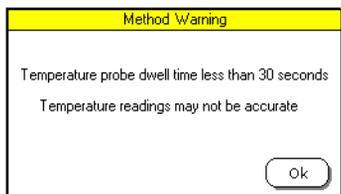
Instrument Calibration Due

This Service Due warning displays when the instrument is due for calibration (“**Menu Screen - Calibration Screen 1**” on page 76). Press **Ok** to clear the warning.



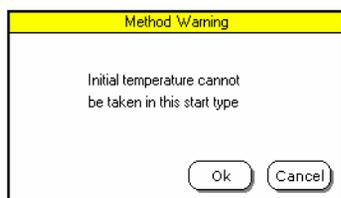
Instrument Preventative Maintenance Due

This Service Due warning displays when the instrument is due for preventative maintenance (“**Preventive Maintenance**” on page 102). Press **Ok** to clear the warning.



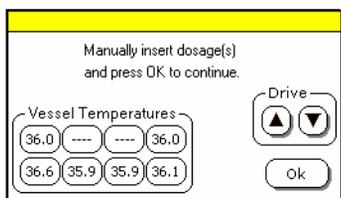
Temperature Probe Dwell Time Warning

This Method Warning screen displays when an Auto Sampling Manifold is installed and enabled (“**Method Properties Screen 3**” on page 62). If the sum of the Manifold Down Duration and Manifold Lead Time (“**Method Properties Screen 3**” on page 62) are less than 30 seconds, this screen displays. Press **Ok** to clear the warning.



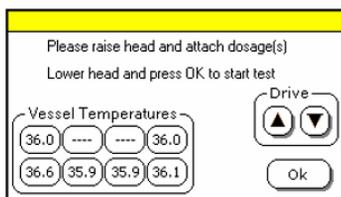
Invalid Initial Temperature Warning

This Method Warning screen displays only when Apparatus 1 (basket) or Apparatus 6 (rotating cylinder) is installed and a Bath Temperature Start is selected from the Start Options screen (“**Method - Start Options**” on page 44). Press **Ok** to clear the warning and start the method without the initial temperature being taken. Press **Cancel** to abort the run and return to the Main screen.

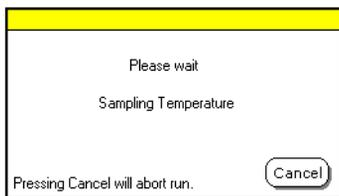
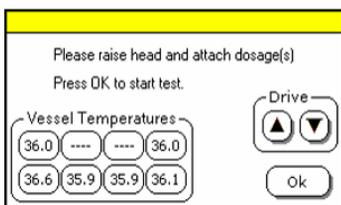


Load Dosage Prompt

This first warning displays when Apparatus 2 (paddle) or Apparatus 5 (paddle over disk) is being used. It indicates that it is time to lower the samples into the vessels. Press **Ok** to clear the dialog and continue the run. Drive ▲ ▼ displays only on units equipped with an automated drive unit.

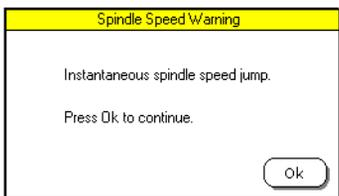


The next two warnings are applicable to systems configured as Apparatus 1 (basket) or Apparatus 6 (rotating cylinder). They indicate that it is time to attach the samples to the shafts. A manual drive unit displays the second warning and an apparatus with an automated drive unit displays the third warning. Press **Ok** to clear the dialog and continue the run. Drive ▲ ▼ displays only on units equipped with an automated drive unit.



Temperature Sampling Notification

This warning displays only when an Auto Sampling Manifold is installed. The warning displays while the temperature probes are immersed in the vessel media and clears itself once the temperature has been read. Press **Cancel** to abort the method and return to the Main screen.

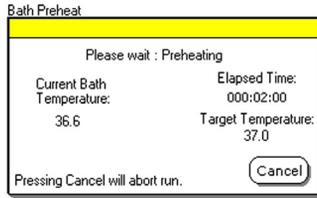


Instantaneous Spindle Speed Jump

If the spindle speed varies by more than 2% over a 40-msec interval, the Spindle Speed Warning screen displays.

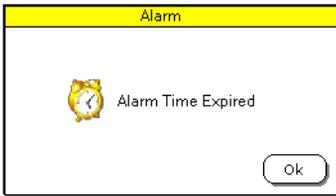
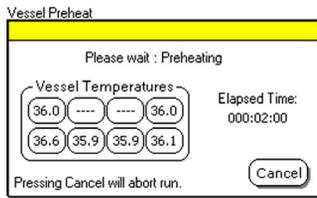
Caution (Stop): The most common cause of this error is a shaft being held while performing diagnostics.

Press **Ok** to return to the Diagnostics screen.



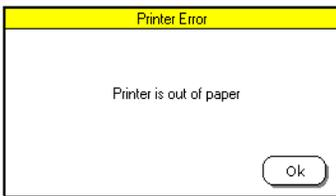
Preheating Notification

This warning displays during preheating for Vessel or Bath selected start options. The warning changes to reflect the current configuration of the system. Current bath or vessel temperatures are displayed during preheating. Press **Cancel** to abort the run and return to the Main screen.



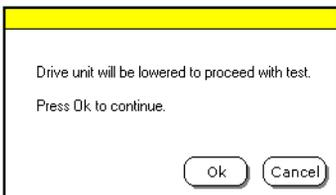
Alarm Time Expired

Displays when the Timer Alarm has expired ("**Main Screen - Alarms**" on page 57). Press **Ok** to clear the warning.



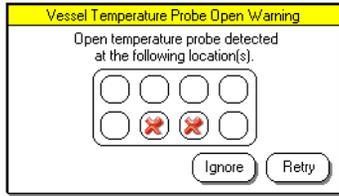
Out of Paper Notification

This warning is applicable to systems with a printer installed. This warning does not interrupt a running method. If the printer runs out of paper during a run ("**Printer Maintenance**" on page 112), the results may be printed out by selecting **Menu > Reports > Last Run Results** ("**Menu Screen - Reports**" on page 67). Press **Ok** to clear the warning.



Automated Drive Unit Movement Warning

Applies to systems with automated drive units. It indicates that you must move the drive unit to its lowest position to continue a run. Press **Ok** to lower the drive unit to its lowest position. Press **Cancel** to abort the current method and return to the Main screen.

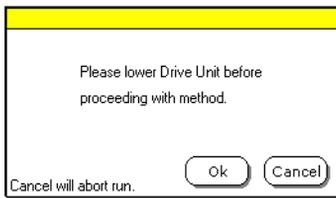


Temperature Probe Open Warning

Indicates that one or more temperature probes have failed. A red X is placed in the appropriate space indicating which probe(s) have failed.

Note: The unit may require service. Refer to the *Agilent 708-DS Service Manual* for troubleshooting.

Press **Retry** to re-test the probes. Press **Ignore** to clear the warning.



Lower Drive Unit Notification

Displays on systems with manual drive units. It indicates that the drive unit must be fully lowered before any testing may proceed. Press **Ok** to retest for the fully lowered position. If the drive unit is not fully lowered, the apparatus beeps and the warning remains on screen. Once the drive unit is in the fully lowered position, press **Ok** to continue the run. Press **Cancel** to abort the current run.



Raise Drive Unit Notification

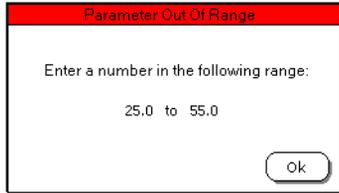
Displays on systems with manual drive units. It indicates when it is permissible to raise the drive unit. Press **Ok** to continue a run. Press **Cancel** to abort the current run.



Delete User

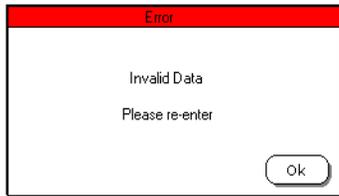
The Delete User screen appears when the Delete User button is pressed on the User Configuration screen. Only an Administrator has the permission to delete 708-DS users.

Error Screens (Red)



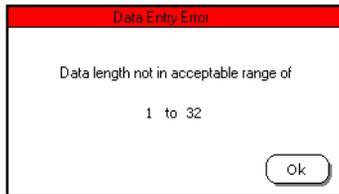
Parameter Out of Range

The Parameter Out of Range error displays when a parameter has been entered that is outside the valid range. The range minimum and range maximum are displayed. This provides an acceptable range and resolution of the data. Press **Ok** to clear the error.



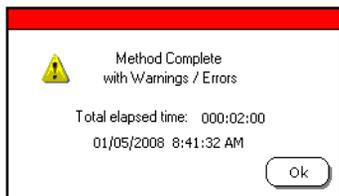
Invalid Data Error

The Invalid Data Error screen displays when unexpected data is encountered. Press **Ok** to clear the error.



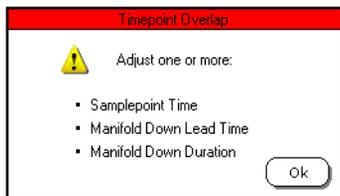
String Length Exceeded

The Data Entry Error displays when a parameter has exceeded the maximum number of characters for that field. The minimum and maximum number of characters are displayed. Press **Ok** to clear the dialog.



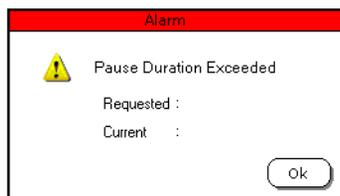
Method Complete with Errors

This error displays after a method has completed and warnings or errors have occurred. The warnings or errors are presented through dialog screens. If a printer is installed, the warning or errors are printed. Press **Ok** to clear the dialog.



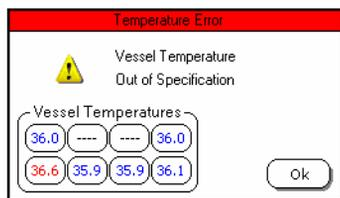
Timepoint Overlap Error

The Timepoint Overlap error screen displays when the timepoint start times entered do not provide sufficient time for all operations to perform before the next timepoint is due. This error typically occurs when the manifold lead time plus the manifold down time exceeds the timepoint interval. See **“Method Properties Screen 4”** on page 64 to correct the overlap. Press **Ok** to clear the error.



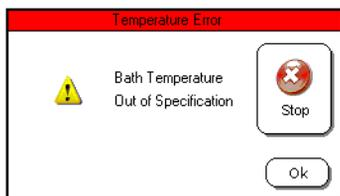
Pause Duration Exceeded

The Alarm error screen displays when the pause duration alarm time has been exceeded (**“Menu Screen - Alarms”** on page 70). Press **Ok** to clear the dialog.



Vessel Temperature Error

Displays when a vessel temperature has exceeded the allowable tolerance range (**“Menu Screen - Instrument Settings Screen 2”** on page 72). The vessel temperatures are displayed in blue if they are within tolerance. They are red if they are outside the allowed tolerance range. Press **Ok** to continue the run.



Bath Temperature Error

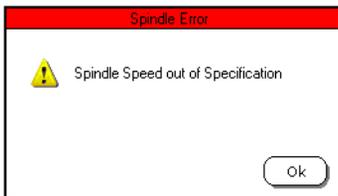
The Temperature Error screen displays when the bath temperature has exceeded the allowable tolerance range (**“Menu Screen - Instrument Settings Screen 2”** on page 72). Press Stop to abort the current method and return to the Main screen. Press **Ok** to continue the run.



Bath Heater Error

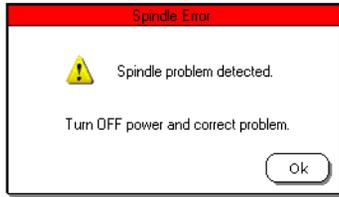
The Bath Error - Heater Disabled screen shows all of the possible errors associated with the bath heater and the bath probe. A red X displays before the current error condition. If a bath heater error is encountered, the current running method is not stopped and you are given the option of correcting the error and continuing the test. Press **Retry** to retest for the error. If the error has been corrected, the dialog clears and you are returned to the Main screen. Press **Ignore** to ignore the error until the unit is powered on / off again.

Bath Heater Error	Action
Heater Probe Open or No Comm.	<ul style="list-style-type: none"> Ensure the bath heater cable is connected. Check cable.
Pump Failed or Heater Off	<ul style="list-style-type: none"> Contact Agilent Customer Care ("Obtaining Warranty and Other Services" on page 116).
Bath Temperature 2 °C Over Setpoint	<ul style="list-style-type: none"> Check for a faulty probe. Contact Agilent Customer Care ("Obtaining Warranty and Other Services" on page 116).
Bath Overtemperature	<ul style="list-style-type: none"> Turn off the bath heater. Contact Agilent Customer Care ("Obtaining Warranty and Other Services" on page 116).
Bath Probe Open	<ul style="list-style-type: none"> Ensure the bath probe is plugged in. Replace the bath probe.
Bath Probe Short	<ul style="list-style-type: none"> Replace the bath probe. Contact Agilent Customer Care ("Obtaining Warranty and Other Services" on page 116).
Heater Probe Short	<ul style="list-style-type: none"> Contact Agilent Customer Care ("Obtaining Warranty and Other Services" on page 116).



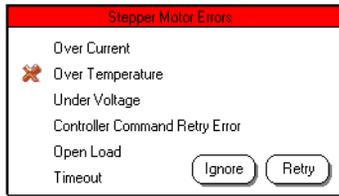
Spindle Speed Error

This Spindle Error screen displays when the spindle speed has exceeded the allowable tolerance range ("**Menu Screen - Instrument Settings Screen 2**" on page 72). Press **Ok** to resume the test and send a message to the printer, if installed.



Fatal Spindle Error

- This Spindle Error screen displays when a serious error has been detected with the spindle motor. For the safety of the equipment and the operator, remove power before correcting the problem. Errors may include jammed paddles, obstructions in the path of the drive belt, and severe deterioration of the spindle motor.
- Contact Agilent Customer Care (“**Obtaining Warranty and Other Services**” on page 116).



Autosampling Manifold Errors

The Stepper Motor Errors screen shows all of the possible errors associated with the manifold stepper motor. A red X displays before the current error condition. If a stepper motor error is encountered, the current run is stopped. You can press **Ignore** to disregard the error, but it must be corrected before a run may be restarted. Press **Retry** to re-test the stepper motor for errors.

Autosampling Manifold Error

Action

Over Current

Over Temperature

Under Voltage

Controller Command Retry Error

Open Load

Timeout

- Contact Agilent Customer Care (“**Obtaining Warranty and Other Services**” on page 116).



5

Maintenance and Troubleshooting

Preventive Maintenance	102
Sample Line / Paddle / Basket Care	103
Vessels and Evaporation Covers	107
Water Bath and Temperature Probes	107
Manifold Maintenance	110
Accessing the Top Cover	111
Printer Maintenance	112

Preventive Maintenance

Preventive maintenance intervals may vary depending on the frequency of instrument usage.

Weekly Maintenance

Inspect the water bath and tubing for algae or other growth. If algae is present, change the water and add algaecide. Ensure the algaecide is compatible with PETG and vinyl plastics.

Monthly Maintenance

If necessary, drain the water from the water bath and clean it thoroughly. Refill the bath and add algaecide.

CAUTION

Ammonia, bleach, alcohol, and other solvents can cause deterioration of the plastic. Use only cleansers approved for plastic materials. These can be found in the catalogs of most scientific supply houses. Additionally, do not use abrasive cleansers that may scratch the plastic.

Bi-Annual Maintenance

- 1 Remove the top cover.
- 2 Inspect the spindle drive belt for wear, fraying, and tension.

NOTE

A small dusting of belt material may be found on the motor plate. This is normal and not a cause for concern.

- 3 Check that the spring-loaded belt tensioner is maintaining pressure on the belt and is not loose.
- 4 Following ESD guidelines, inspect all printed circuit board assemblies (PCBA) for corrosion and debris. If debris is found, use pressurized inert gas to expel it from the board. Also, ensure that the PCBAs are securely attached.
- 5 Following ESD guidelines, ensure all connectors are securely fastened to each PCBA and to the power supply, switch, sensors, or motors. Also inspect the connectors and cabling for damage.
- 6 Lubricate the linear bearing upright supports with 3-in-1 oil.

Sample Line / Paddle / Basket Care

Shafts

- When using with corrosive materials such as hydrochloric acid or media containing salts, be sure to rinse them thoroughly with deionized water immediately after each use, and dry thoroughly with a soft towel or cloth.
- Do not clean with abrasive cleansers or cloths. Use deionized water whenever possible. If you must use a cleanser or solvent, be sure that it is as mild as possible, non-abrasive, and fully compatible with fluorocarbons and stainless steel before use. If in doubt, call the service department for advice before proceeding.
- We recommend that you do not use a laboratory dishwasher. Clean paddles and basket shafts only by hand. The high temperatures to which your items would be subject in a dishwasher may damage the fluorocarbon coating.
- Be sure to handle with care. If you must clean or handle the shafts while they are still mounted on the instrument, use minimal pressure on the shaft to prevent them from bending. When secured in the drive unit of the instrument, just a little bit of pressure exerted on the shaft—especially near the blade or basket—can easily bend the shaft and cause significant wobble.
- Use care when removing vessels from the apparatus while the paddles or basket shafts are installed so that you do not bump them.
- When attaching or removing baskets, do not bend the clips excessively.

- Please store paddles and basket shafts properly between uses. Do not simply place these items in a drawer. They will be subject to nicks, chips, and scratches as they bump against each other. Place them back into the original styrofoam shipping container or other appropriate container between uses. This will prevent them from coming into contact with each other or anything else in the storage area.

NOTE

Place protective caps on dry and clean threads when storing spin on/off shafts.

Baskets

- When using with corrosive materials such as hydrochloric acid or media containing salts, be sure to rinse them thoroughly with deionized water immediately after each use, and dry thoroughly with a soft towel or cloth.
- Please do not clean baskets with abrasive cleansers or cloths, especially if they're gold or PTFE coated (non-stick). Mesh openings on baskets could enlarge, which could have an effect on results. Use deionized water whenever possible. If you must use a cleanser or solvent, be sure that it is as mild as possible, non-abrasive, and fully compatible with fluorocarbons and stainless steel before use. If in doubt, contact the service department for advice before proceeding.
- We recommend that you do not use a laboratory dishwasher. Clean baskets only by hand. The high temperatures to which your baskets would be subject in a dishwasher may damage the fluorocarbon coating.
- Use caution when handling baskets. It is important that they retain their cylindrical shape, so take care not to kink or bend the mesh. Check frequently to ensure that the mesh is completely open and that there are no rips or tears.
- Please store baskets properly between uses. Do not simply place these baskets in a drawer. They will be subject to nicks, chips, and scratches as they bump against each other and they may get bent out of shape. Place them back into the original shipping container or other appropriate container between uses. This will prevent them from coming into contact with each other or anything else in the storage area.

Rinsing Sample Lines

If the Agilent 708-DS is configured with a sampling manifold, the sample lines require a routine rinsing after each test. A rinse tubing kit (17-1341) or rinse tray is provided to conveniently flush the sample tubing.

Rinse Tubing Kit

- 1 Attach one end of the lines of the rinse tubing kit to the sample cannulas and place the other in a container of an appropriate rinsing solution.
- 2 Activate the pumping mechanism to properly flush the tubing lines.

Rinse Tray

- 1 Raise the drive unit of the 708-DS to its maximum height.
- 2 Push each shaft up to provide clearance for the rinse tray.

NOTE

Do not rotate the shafts with the shafts pushed up. This could cause damage to the cannulas / probes on the sampling manifold.

- 3 Fill the rinse tray with the desired cleaning solution.
- 4 Position the rinse tray on the vessel plate so that each sample cannula is above the tray.
- 5 Lower the sampling manifold.
- 6 Carefully lower the drive unit of the 708-DS until each sample cannula is submerged in the cleaning solution of the rinse tray.
- 7 Activate the pump configured with the system until the tubing is sufficiently cleansed.
- 8 Raise the manifold and drive unit.
- 9 Remove the rinse tray and properly dispose of the rinsing solution.
- 10 Lower the shafts to their appropriate location.

Rinse Cups (Optional)

If the Agilent 708-DS is configured with a sampling manifold, the sample lines require a routine rinsing after each test. Easy-to-attach rinse cups are available to make the flushing of the tubing simple and efficient.

- 1 Fill the rinse cups with the cleaning solution desired and attach a cup to each shaft.



Figure 66. Rinse Cup

- 2 The sample cannulas must be sufficiently immersed in the solution. The return cannulas can either be contained in the cup or allowed to drip into the vessel. The bottom of the cup rests on either the paddle blade or the basket clip attachment.
- 3 Activate the pump configured with the system until the tubing loop is sufficiently cleansed.
- 4 Finally, detach the rinse cups, place them in their transport tray, dispose of the waste, and rinse the cups to be ready for their next use.

Vessels and Evaporation Covers

- 1 Carefully wash the dissolution vessels after each use.

Water Bath and Temperature Probes

Water Bath Care

The water bath supplied with the 708-DS is designed to be maintenance free except for occasional cleaning.

CAUTION

Ammonia and bleach can cause deterioration of the plastic. Use only cleansers approved for plastic materials. These can be found in the catalogs of most scientific supply houses. Additionally, do not use abrasive cleansers that may scratch the plastic.

If algacide or a clear bath product is used, ensure that it is compatible with PETG and vinyl. The flow path in the heater / circulator is primarily stainless steel and should tolerate most clear bath formulations. Check with the product manufacturer for compatibility.

Complete these steps if any of the water bath fittings are leaking:

- 1 Turn off the heater / circulator and drain the water bath completely.
- 2 If the inlet to the bath is leaking, remove the leaking right-top bath bulkhead fitting and unscrew the elbow fitting.
- 3 If the outlet to the bath is leaking, remove the tubing and unscrew the T-fitting from the left-bottom side of bath.
- 4 Inspect the bulkhead fitting gaskets for damage and replace them if necessary.

- 5 Remove old PTFE tape from all male fittings. Inspect threads for damage and replace the elbow of the T-fitting if necessary.
- 6 Apply new PTFE tape to the male fitting threads.
- 7 Reinstall and tighten the bulkhead fitting to the bath and / or screw in the elbow or T-fitting, depending on whether it's an inlet or outlet.
- 8 Fill the bath with water and inspect for leaks. If it's still leaking, contact Technical Service.

Water Bath Temperature Probe Accuracy Test

The temperature probe can be tested using a high-quality ohmmeter capable of at least four-digit resolution and a known temperature bath. Both the multi-meter and the bath temperature must be traceable to a known reference standard such as NIST.

- 1 Place the probe into a water bath with a known temperature and allow several minutes for the probe to equilibrate.
- 2 The resistance is measured by attaching the leads of the multi-meter to the top and shank of the ¼-inch phone plug.
- 3 Refer to the table on Page [page 109](#) to find the resistance value of the probe at the bath temperature. Temperatures between the values listed may be interpolated.
- 4 Probes are interchangeable and manufactured with a tolerance of +/- 0.1 °C. Probes found to be outside of the tolerance of +/- 0.1 °C should be replaced.

Temp. Centigrade	Resistance Ohms	Temp. Centigrade	Resistance Ohms	Temp. Centigrade	Resistance Ohms
-5.0	9530	24.0	2354	53.0	724.50
-4.0	9046	25.0	2252	54.0	697.90
-3.0	8586	26.0	2156	55.0	672.50
-2.0	8151	27.0	2064	56.0	648.10
-1.0	7741	28.0	1977	57.0	624.80
0.0	7355	29.0	1894	58.0	602.40
1.0	6989	30.0	1815	59.0	580.90
2.0	6644	31.0	1739	60.0	560.30
3.0	6319	32.0	1667	61.0	540.50
4.0	6011	33.0	1599	62.0	521.50
5.0	5719	34.0	1533	63.0	503.30
6.0	5444	35.0	1471	64.0	485.80
7.0	5183	36.0	1412	65.0	469.00
8.0	4937	37.0	1355	66.0	452.90
9.0	4703	38.0	1301	67.0	437.40
10.0	4482	39.0	1249	68.0	422.50
11.0	4273	40.0	1200	69.0	408.20
12.0	4074	41.0	1152	70.0	394.50
13.0	3886	42.0	1107	71.0	381.20
14.0	3708	43.0	1064	72.0	368.50
15.0	3539	44.0	1023	73.0	356.20
16.0	3378	45.0	983.80	74.0	344.50
17.0	3226	46.0	946.20	75.0	333.10
18.0	3081	47.0	910.20	76.0	322.30
19.0	2944	48.0	875.80	77.0	311.80
20.0	2814	49.0	842.80	78.0	301.70
21.0	2690	50.0	811.30	79.0	292.00
22.0	2572	51.0	781.10	80.0	282.70
23.0	2460	52.0	752.20		

Figure 67. Resistance Value of Temperature Probe

Removing or Adjusting Cannulas / Probes

The manifold of the apparatus may contain up to two cannulas and one probe per vessel position depending on the options configured. To adjust or remove these items:

- 1 Loosen the thumbscrew on the individual position of the manifold and modify the location of the cannula / probe accordingly.

Manifold Maintenance

Removing the Manifold

The manifold of the Agilent 708-DS is installed and removed with the touch of a button. The initial installation procedure is described in **“Sample Manifold Installation”** on page 27.

For the purposes of cleaning, maintenance, or just to create additional space for manual testing, the manifold may be ejected from the drive unit. To remove the manifold, follow these steps:

- 1 Lift the drive unit to its fully raised position.
- 2 Disconnect the two thermistor RS232 cables (if AutoTemp is installed).
- 3 Remove the sample tubing from the black fasteners.
- 4 Remove the evaporation covers.
- 5 Balance the manifold with one hand and press **Eject** to fully lower the manifold until it is free from the drive unit. Catch the manifold as it is released.

The Manifold Control screen can be accessed by pressing **Menu > Diagnostics**.

Accessing the Top Cover

Removing Top Cover

- 1 Ensure the Lift Drive is in the full, down operational position. Turn off the power and remove the power cord.
- 2 Lower all paddle / basket shafts or remove them from the machine.
- 3 Remove the top screws from the front and back of the cover. Remove the left- and right-side screws from bottom sides of cover.
- 4 If a printer is installed in the top cover, lift the cover up approximately six inches and remove the two cable connectors from the printer inside.
- 5 Remove the cover away from the instrument.

Replacing the Top Cover

- 1 If a printer is installed in the top cover, reconnect the cable connectors to the printer inside the cover.
- 2 Press the cover down into place, aligning the front holes with the respective dosage delivery module (DDM) openings, and slipping over the left-hand power switch.
- 3 Re-fasten the cover screws.

Printer Maintenance

Removing a Thermal Paper Roll

NOTE

A red line on the side of the paper roll indicates the paper supply is almost exhausted. Replacement is recommended when the red line is visible. If the paper roll is not changed and the paper supply is exhausted, the 708-DS gets an out-of-paper signal from the printer, and displays an error on the screen.

- 1 Place fingers on each side of the printer panel and pull forward to expose the paper chamber and the paper roll, or empty core inside.



Figure 68. Accessing the Printer Handle

- 2 Use your fingers to remove the small roll or core.

NOTE

The empty core is disposable. Each new roll includes a cardboard core.

Inserting a Thermal Paper Roll

NOTE

Only a 58 mm wide x 13.65 meter (maximum) long thermal printer paper roll can be used in the printer. For an FDA environment, it must also meet the ten-year retention requirement. Ordering replacement paper from Agilent is recommended.

- 1 With the paper chamber exposed, insert the paper roll so the leading edge of the paper will feed over the top of the printer panel.



Figure 69. Inserting a Thermal New Paper Roll

- 2 Keeping the paper centered, close the printer panel and snap it into place.



Figure 70. Closing the Printer Chamber

- 3 To verify that the paper is not skewed or jammed, press the button on the top-left of the printer panel to feed some paper. If paper does not feed, re-open the panel and re-center the paper.



Figure 71. Feeding Paper

Thermal Printer Test

In the second Diagnostics screen, press **Print Test Sheet** to print sample text and verify the printer is functioning properly.

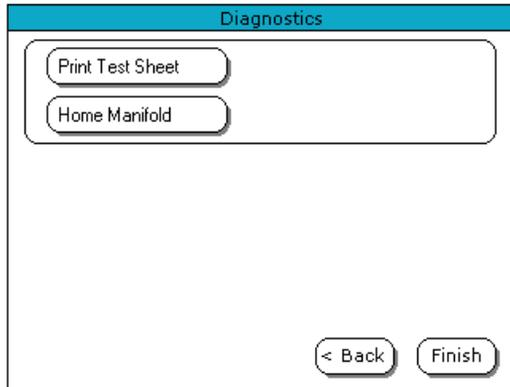


Figure 72. Printer Test

Obtaining Warranty and Other Services

To place a service order (warranty or other services), please contact your local Customer Care Center. Contact information can be found at www.agilent.com under your country using the Contact Us link. Place your service request using the displayed phone number or E-mail address.

In This Book

- Chapter 1 Safety
- Chapter 2 Introduction
- Chapter 3 Setup
- Chapter 4 Operation
- Chapter 5 Maintenance and Troubleshooting

www.agilent.com

© Agilent Technologies, Inc. 2020

70-9058 Rev H February 2020

