

750
Heater / Circulator

Operator's Manual



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2011

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

Manual Part Number

70-9063

Edition

Rev A, March 2011

Printed in USA

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.

Contents

| | | |
|----------|---|-----------|
| | Figures | 5 |
| | Tables | 7 |
| 1 | Safety | 9 |
| | Electrical Hazards | 9 |
| | Warning | 11 |
| | Caution | 11 |
| | Note | 11 |
| | Information Symbols | 12 |
| | General | 13 |
| | WEEE Directive | 14 |
| 2 | Introduction | 15 |
| | Conventions Used in this Manual | 17 |
| 3 | Setting Up the 750 | 19 |
| | Unpacking Your 750 | 20 |
| | Electrical Connections | 21 |
| | Interfacing to the Dissolution Apparatus | 22 |
| | Liquid Connections—Single Apparatus Systems | 23 |
| | Liquid Connections—Dual Apparatus Systems | 25 |
| | Hidden Key Functions | 27 |
| 4 | Operating the 750 | 29 |
| | Initial Power Up | 30 |

Contents

| | |
|---|-----------|
| Priming the System | 31 |
| Operation Under Apparatus 1 / 2 Control | 32 |
| Set Water Bath Temperature | 32 |
| Set Time Delay Start | 32 |
| Operation Under Apparatus 3 / Apparatus 7 Control | 33 |
| Set Water Bath Temperature | 33 |
| Set Time Delay Start | 33 |
| Operation Under Local Control | 34 |
| Over Temp | 35 |
| Optional Report Center Printer | 36 |
| Setting Date and Time | 37 |
| Setting Automatic Printout Interval | 37 |
| RS232 Command Set | 38 |
| RS232 Connections | 39 |
| 5 Maintenance | 41 |
| Periodic Maintenance | 42 |
| Monthly | 42 |
| Fuse Replacement | 43 |
| 6 Service and Warranty | 45 |
| Service and Warranty Information | 46 |
| Exclusions and Limitations | 46 |
| Obtaining Warranty Service | 46 |
| Warranty Limitations | 47 |
| Exclusive Remedies | 47 |
| Index | 49 |

Figures

- Figure 1. 750 Heater / Circulator 16
- Figure 2. 750 Rear Panel 21
- Figure 3. Single Water Bath Connections 23
- Figure 4. Dual Water Bath Tubing Connections 25
- Figure 5. 750 Keypad 34
- Figure 6. Report Center Printer 36
- Figure 7. DB-9 Connector for RS232 39

Figures

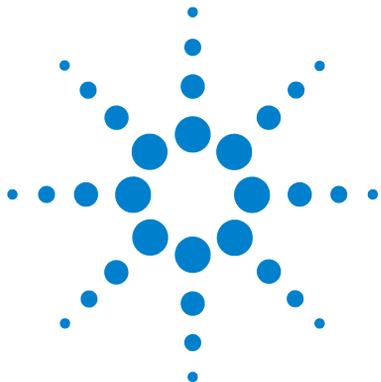
This page was intentionally left blank, except for this message.

Tables

| | |
|-------------------------------|----|
| Table 1. Hidden Key Functions | 27 |
| Table 2. RS232 Command Set | 38 |

Tables

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



1 Safety

Electrical Hazards 9
WEEE Directive 14

The 750 has been carefully designed 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 a 750 involves the use of aqueous liquids. 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 750 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 8020 to determine which parts are operator-accessible.



1 Safety

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 consensus-based standards of safety to provide assurance, required by OSHA, that 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

General

CE Compliant Products

The 750 has been designed to comply with the requirements of the Electro-magnetic Compatibility (EMC) Directive and the Low Voltage Directive (LVD) of the EU.

Agilent, Inc. has confirmed that each product complies with the relevant directives by testing a prototype against the prescribed European Norm (EN) standards.

Proof that a product complies with the directives is indicated by:

- the CE marking appearing on the rear of the product.
- the documentation package that accompanies the product containing a copy of the declaration of conformity. This declaration is the legal declaration by Agilent, Inc. that the product complies with the directives and also shows the EN standards to which the product was tested to demonstrate compliance. The declaration of conformity is signed by the representative of the manufacturing plant.

cTUVus - U.S. and Canadian Product Approvals

The 750 has been designed to comply with North American safety requirements.

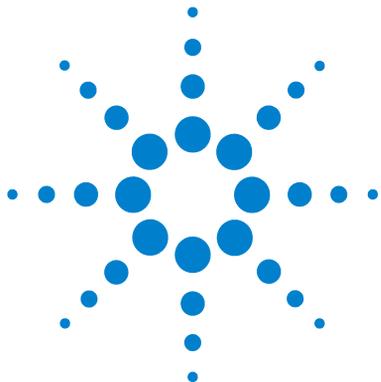
This product has been tested and certified for the North American market by TUV Rheinland of North America, Inc. The TUVus mark signifies that this product has been tested to U.S. standards and certified for the U.S. market. The cTUV mark signifies that this product has been tested to Canadian standards and certified for the Canadian market. When the two marks are coupled, the cTUVus mark signifies that this product has been tested to standards and certified for both markets.

WEEE Directive

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 wheellie bin” WEEE symbol shown on page 12 in accordance with European Standard EN 50419.

This symbol on the product or on its packaging indicates that this product must not be disposed of as unsorted municipal waste. The separate collection and recycling of your waste equipment at the time of disposal will help to conserve natural resources and ensure that it is recycled in a manner that protects human health and the environment.

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



2 Introduction

Conventions Used in this Manual 17

The 750 Heater / Circulator has been designed for precise temperature control of your dissolution water bath with a range from ambient to 55 °C and a precision of ± 0.1 °C.

The operating parameter values can be set and monitored either from the front panel of the instrument or with the keypad of your dissolution apparatus. An interface cable interconnects the units enabling them to communicate with each other. The temperature is shown on a large, easy-to-read digital display, and the instrument status is continuously indicated on individual LEDs.

For the highest accuracy, the 750 can be used with an optional plug-in remote temperature probe for direct monitoring of the water bath temperature. This probe is supplied with all VK 7000 / 7010 Dissolution Apparatus as a standard accessory. As an alternative, the temperature probe inside the internal tank is used for temperature control. For easy automation, all functions can be controlled and monitored via RS232 with both DB-9 and RJ-11 connectors provided on the rear panel.

WARNING

The equipment 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.



The optional Report Center Printer gives you a printout of the system status at user-selectable time intervals for easy documentation of operations. This feature is designed to help you comply with regulatory requirements.



Figure 1 750 Heater / Circulator

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

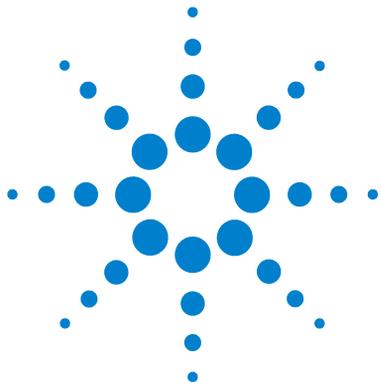
- Items you are asked to press are in bold. For example, “press **1** on the keypad”.
- Key sequences you are asked to press appear like this: **CLEAR > 7**.

NOTE

Remember to return the warranty card. Completing and returning the card ensures your right to protection under the terms and conditions of your warranty. It also enables us to better assist you in the event of any problems. Additionally, it guarantees you will be informed of any issues arising concerning your equipment, such as upgrades, retrofits, or regulatory changes.

2 Introduction

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



3 Setting Up the 750

| | |
|---|----|
| Unpacking Your 750 | 20 |
| Electrical Connections | 21 |
| Interfacing to the Dissolution Apparatus | 22 |
| Liquid Connections—Single Apparatus Systems | 23 |
| Liquid Connections—Dual Apparatus Systems | 25 |
| Hidden Key Functions | 27 |



Unpacking Your 750

Complete the following steps to unpack your 750:

- 1 Carefully remove all items from the shipping carton.
- 2 Check all items for damage during shipping. If any damage to the instrument is evident, contact both the carrier who delivered the instruments to you and Agilent, Inc. Though claims for damage should be filed with the carrier, we will be glad to help you in filing a claim and in getting your system up and running as quickly as possible.
- 3 Check the shipping carton for any items which may have come loose during shipping before discarding or storing the packaging.
- 4 Place the unit on a clear, dry, and level section of the benchtop as close to the dissolution apparatus as possible. At least eight inches (20 cm) of unobstructed space should be available behind the unit for easy access to the rear panel power and liquid connections. As with any electronic apparatus, the area around the instrument must be kept clean and dry.

WARNING

The electrical connection at the back of the equipment is the primary disconnect for the instrument. The heater / circulator should be positioned to allow accessibility to the power cords for easy disconnection.

A properly grounded, GFC recommended AC power receptacle rated at 20 amps or higher should be available within six feet (two meters) of the unit.

NOTE

If you are interfacing your 750 to the VK 7000 / 7010 Dissolution Apparatus, you can place it lengthwise in the space under the rear of the vessel table behind the molded water bath. Situate the heater / circulator so that you have access to the AC power switch (the keypad is not accessible). Should you need to move the heater / circulator, it is necessary to lift the entire dissolution apparatus. Therefore, this option is recommended only if bench space is at a premium.

Electrical Connections

WARNING

Ensure the equipment is configured for the voltage supplied.

Complete the following steps to make electrical connections:

- 1 Ensure the power switch on the rear panel AC power connector is in the OFF position. See Figure 2, “750 Rear Panel,” below.



Figure 2 750 Rear Panel

- 2 Connect the power cord to the AC power connector.

WARNING

Because of the wide variety of AC outlet formats around the world, the power cord you need may not be the one supplied with the unit. It may be necessary for you to supply the power cord needed for your country or area. Be sure that any power cord you use meets all local safety regulations and is rated to carry at least 10 amps.

- 3 Plug the other end of the cord into an outlet of the proper voltage.

Interfacing to the Dissolution Apparatus

If you are using your 750 Heater / Circulator with a VK 7000 / 7010 Dissolution Apparatus or Apparatus 3 / Apparatus 7, you can control the 750 with the dissolution apparatus keypad. To interface the units, complete the following steps:

- 1 Locate the four-pin cable in the 750 accessory kit. Plug one end into the 750 rear panel jack labeled INPUT SIGNAL and the other end into the jack on the dissolution apparatus rear panel labeled HEATER/CIRC.

NOTE

When the 750 is interfaced to the VK 7000 / 7010, the heater / circulator LED displays a series of dashes (---), indicating it is under remote control and the keypad is locked.

- 2 Ensure the remote temperature probe supplied with the dissolution apparatus is plugged into the rear panel jack labeled BATH TEMP. This enables the dissolution apparatus to monitor and display the water bath temperature while controlling the heater / circulator.
- 3 Set the desired water bath temperature via the dissolution apparatus Main Menu option 3 (Set Temperature on the VK 7000 / 7010 or Bath Temperature on the Apparatus 3 / Apparatus 7).

Liquid Connections—Single Apparatus Systems

Use the following figure and the steps below to connect a single dissolution apparatus water bath to the 750:

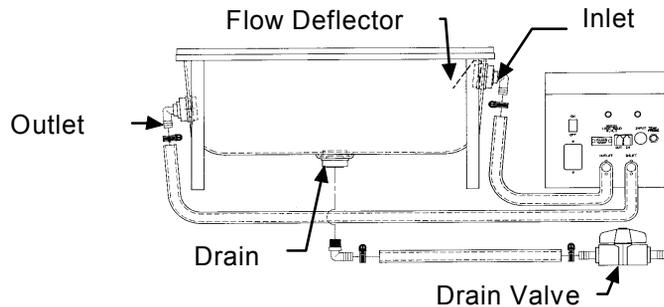


Figure 3 Single Water Bath Connections

- 1 Locate the two pieces of plastic tubing and the four stainless steel tubing clamps included in the 750 tubing kit.
- 2 Slip one end of each piece of tubing over the barbed 90-degree angle adapters in the inlet and outlet bulkhead fittings of the water bath. The inlet fitting is the higher of the two fittings. It has a perforated deflector screen attached on the inside of the water bath (see [Figure 3, “Single Water Bath Connections,”](#) above). Slide the tubing as far up on the adapter as possible.
- 3 Slip a tubing clamp over the free end of each piece of plastic tubing and secure each clamp over the barbed angle adapters. To avoid damage to the tubing, do not overtighten.
- 4 Slip the other pair of tubing clamps over the free ends of each piece of tubing and insert the tubing ends onto the inlet and outlet connectors on the rear of the 750. Ensure the tubing connected to the water bath inlet attaches to the 750 rear panel connector labeled OUTLET and the tubing connected to the water bath outlet attaches to the 750 rear panel connector labeled INLET. Secure the clamps over the inlet and outlet connectors. To avoid damage to the tubing, do not overtighten.
- 5 Double check all connections.

3 Setting Up the 750

- 6 Fill the water bath with purified water to approximately one inch (2.5 cm) above the inlet fitting. The system is now ready for use.

NOTE

Use ultrapure water when possible to minimize scale and mineral buildup. Use algacide to inhibit mold and bacteria 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 750 are primarily stainless steel and should tolerate most clear water bath formulations.

Liquid Connections—Dual Apparatus Systems

When using your 750 with two dissolution apparatus configured as 12-, 14-, or 16-spindle systems, the 750 output is divided between the two water baths, then recombined at the inlet. The water baths are also connected via the drains to ensure the water levels are always equal.

Use the following figure and the steps below to connect a dual dissolution apparatus water bath to the 750:

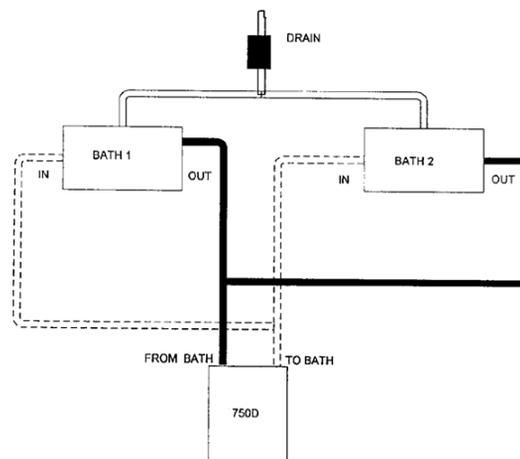


Figure 4 Dual Water Bath Tubing Connections

- 1 Locate the two short lengths of tubing included in the 750 tubing kit.
- 2 Slip one end of each piece of tubing over the inlet and outlet connectors on the 750 rear panel. Slip a tubing clamp over the free end of each piece of plastic tubing and secure each clamp over the connectors. To avoid damage to the tubing, do not overtighten.
- 3 Slip a tubing clamp over the free end of each piece of tubing and insert one of the supplied Y connectors into the end of each. Secure the tubing clamps over the Y connectors. To avoid damage to the tubing or connectors, do not overtighten.

3 Setting Up the 750

- 4 Attach a length of tubing to both branches of the Y connector on the 750 outlet. Slip a tubing clamp over the free end of each piece of tubing and secure the tubing clamp over the Y connector. To avoid damage to the tubing, do not overtighten.
- 5 Slip another tubing clamp over the free end of each piece of tubing and attach the free end of each piece of tubing to the barbed angle adapters on the water bath inlets. The inlet fitting is the higher of the two water bath fittings and has a perforated deflector screen attached on the inside of the water bath. Secure the tubing clamps. To avoid damage to the tubing, do not overtighten.
- 6 Repeat steps 4 - 5 to attach the 750 inlet connectors to the water bath outlets.
- 7 Locate the length of tubing with the drain valve inserted. Slip a tubing clamp over each end and secure the tubing clamps over the drain on the bottom of each water bath. To avoid damage to the tubing, do not overtighten.
- 8 Double check all connections.
- 9 Fill the water baths with purified water to approximately one inch (2.5 cm) above the inlet fitting. The system is now ready for use.

NOTE

Use ultrapure water when possible to minimize scale and mineral buildup. Use algacide to inhibit mold and bacteria 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 750 are primarily stainless steel and should tolerate most clear water bath formulations.

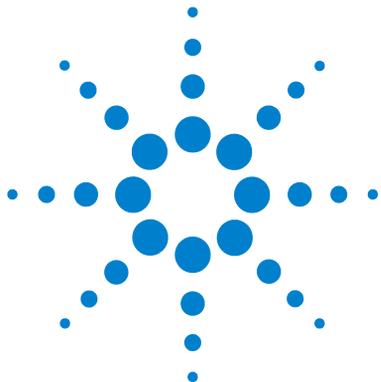
Hidden Key Functions

Table 1 Hidden Key Functions

| Key Sequence | Function |
|---------------------|---|
| CLEAR > 1 | Use this key sequence to toggle on and off the Report Center Printer (if installed). |
| CLEAR > 4 | Use this key sequence to toggle the position of the decimal on the temperature display from one tenth of a degree (0.1 °C) to one one-hundredth of a degree (0.01 °C). NOTE: This function does not require that you press ENTER to complete the task. |
| CLEAR > 7 | Use this key sequence to enter a different communication port identification number. Enter the appropriate identification number and press ENTER . |
| CLEAR > 8 | Use this key sequence to toggle between baud rates of 1200, 2400, 4800, and 9600 BPS. Select the appropriate baud rate by continuing to press CLEAR until the desired selection displays. Press ENTER . |

3 Setting Up the 750

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



4 Operating the 750

| | |
|---|----|
| Initial Power Up | 30 |
| Priming the System | 31 |
| Operation Under Apparatus 1 / 2 Control | 32 |
| Operation Under Apparatus 3 / Apparatus 7 Control | 33 |
| Operation Under Local Control | 34 |
| Over Temp | 35 |
| Optional Report Center Printer | 36 |
| RS232 Command Set | 38 |
| RS232 Connections | 39 |

CAUTION

Before operating the circulator, ensure the water bath is filled and the system is fully primed as described in “[Priming the System](#)” on page 31. Prolonged operation without water damages the pump and heaters.



Initial Power Up

Turn on the heater / circulator. Flow begins almost immediately. The unit runs very quietly with only a low hum. If you hear gurgling noises, it is because air remains in the tubing. The air is usually expelled within approximately 30 seconds. If it persists, stop and prime the system as described in [“Priming the System”](#) on page 31.

Priming the System

The 750 Heater / Circulator has been designed to be self-priming under most conditions. However, it may require some assistance, especially when it is new or has not been used for a period of time, so that the tubing is completely dry.

Complete the following steps to prime the unit:

- 1 Ensure the water bath is adequately filled.
- 2 Turn on the pump and tilt up the rear of the unit. This helps the air inside the unit to flow out with its natural buoyancy. You should see and hear the air being expelled through the outlet tubing.
- 3 If more vigorous intervention is required, hold the unit below the level of the benchtop (if the tubing is long enough to allow you to do so) while tilting up the rear.
- 4 Should the unit fail to prime after these steps, contact the Dissolution Systems Service Department or your local Agilent representative for assistance.

Operation Under Apparatus 1 / 2 Control

Set Water Bath Temperature

- 1 From the system status screen, press **MENU**. The Main Menu displays.
- 2 Select option **3**, Set Temperature. The Select Temperature Set Mode screen displays.
- 3 Select option **1**, Set Bath Temp.
- 4 Enter the desired water bath temperature and press **ENTER**. The newly set temperature displays on the dissolution apparatus. If the actual temperature is lower than the set temperature, the Heater On LED illuminates, indicating the heating element is on.

Set Time Delay Start

- 1 From the system status screen, press **MENU**. The Main Menu displays.
- 2 Select option **3**, Set Temperature. The Select Temperature Set Mode screen displays.
- 3 Select option **3**, Delayed Bath Heating.
- 4 Enter the desired water bath temperature and press **ENTER**.
- 5 Enter the desired start time in 24-hour format and press **ENTER**.
- 6 Enter the desired start date in the appropriate format and press **ENTER**. The system status screen displays.

Operation Under Apparatus 3 / Apparatus 7 Control

Set Water Bath Temperature

- 1 From the Ready screen, press **MENU**. The Main Menu displays.
- 2 Select option **3**, Bath Temperature. The Bath Temperature Control/Report screen displays.
- 3 Select option **1**, Set Bath Temp.
- 4 Enter the desired water bath temperature in xx.x format and press **ENTER**. The Main Menu displays. If the actual temperature is lower than the set temperature, the Heater On LED illuminates, indicating the heating element is on.

Set Time Delay Start

- 1 From the Ready screen, press **MENU**. The Main Menu displays.
- 2 Select option **3**, Bath Temperature. The Bath Temperature Control/Report screen displays.
- 3 Select option **2**, Delay Heating.
- 4 Enter the desired start time in 24-hour format and press **ENTER**.
- 5 Enter the desired start date in the appropriate format and press **ENTER**. The top line of the display continues to flash until the function is cancelled or until the set date and time are reached.
 - To cancel the delayed heating function, press **ESC**. The Ready screen displays.

Operation Under Local Control

If you do not interface your 750 to an Apparatus 1 / 2 or Apparatus 3 / 7, operational parameters are set via the 750 keypad as shown below.

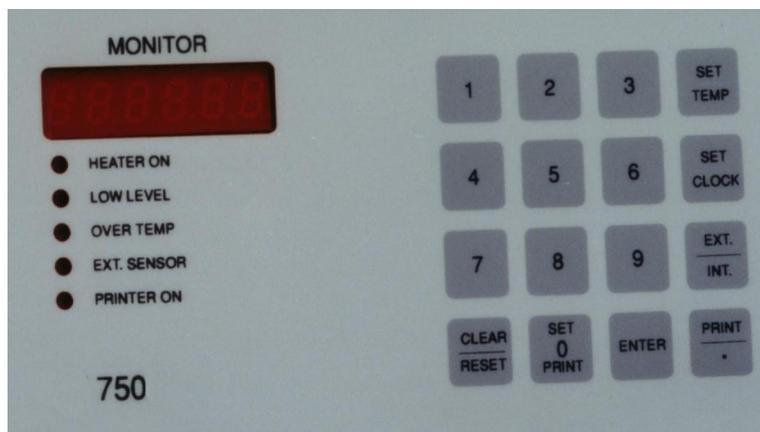


Figure 5 750 Keypad

Complete the following steps to set the operational parameters using the 750 keypad:

- 1 If you have a remote temperature probe plugged into the 750 rear panel, check the Ext. Sensor LED on the front panel. If it is illuminated, the unit is using the remote probe to monitor the actual water bath temperature. If it is not illuminated, the internal sensor is being used. Press **EXT./INT.** to toggle between the internal and external sensors.
- 2 Set the desired water bath temperature by pressing **SET TEMP.** The current setting flashes. Press **ENTER** to accept or key in the desired temperature in xx.x format. Press **ENTER.** The Heater On LED illuminates if the set temperature is higher than the actual temperature and stays on until the water reaches the desired temperature. It flashes intermittently as the heater approaches the set point. The digital display shows the actual temperature of the circulating water on a continuous basis.

Over Temp

The Over Temp LED illuminates to indicate that the heater has exceeded its 55 °C maximum temperature. This rare event may occur if the water level drops suddenly or if a temperature sensor malfunctions. Shut down power to the instrument and troubleshoot the system. If the source of the problem cannot be located, contact the Dissolution Systems Service Department or your local Agilent representative for assistance.

Optional Report Center Printer

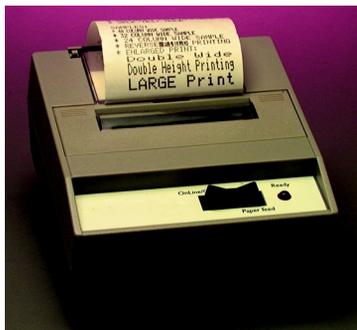


Figure 6 Report Center Printer

NOTE

Use the hidden key function CLEAR > 1 to enable the Report Center Printer prior to programming it for use.

The 750 Heater / Circulator is available with an optional Report Center Printer for use as a standalone machine. It provides hard-copy documentation of the unit's activity, printing the elapsed time, time-of-day, and the actual temperature at user-selectable time intervals.

NOTE

The PRINT key toggles on and off the Report Center Printer. Its status is indicated by the Printer On LED on the front panel.

Setting Date and Time

If applicable, complete the following steps to set the correct date and time:

- 1 Press **SET CLOCK**. The current date flashes in mm/dd/yy format.
- 2 Press **ENTER** to accept the value displayed or enter the correct date and press **ENTER**. The current time flashes.
- 3 Press **ENTER** to accept the value displayed or enter the correct time in 24-hour format (for example, 0900 for 9:00 AM or 1330 for 1:30 PM) and press **ENTER**.

Setting Automatic Printout Interval

If applicable, complete the following steps to set the automatic printout interval:

- 1 Press **SET PRINT**. The current setting flashes in minutes.
- 2 Press **ENTER** to accept the displayed value or enter the new time interval up to 999 minutes and press **ENTER**.

RS232 Command Set

The following RS232 codes are available for external communication and/or control of your 750 Heater / Circulator via the rear panel RS232 serial ports:

Table 2 RS232 Command Set

| Function | Mode | Command | Response | Variables |
|-----------------|-------|------------|------------|---|
| Time | Read | 0123000000 | 012300hhmm | hh = hours mm = minutes |
| | Write | 013300hhmm | Same | |
| Date | Read | 0123100000 | 01231mmddy | mm=month dd = day y = year (last digit) |
| | Write | 01331mmddy | Same | |
| Set Bath Temp. | Read | 0122000000 | 01220fttt | tttt = ttt.t degrees f = 0 Celsius f = 2 Fahrenheit |
| | Write | 01320fttt | Same | |
| Actual Temp. | Read | 0122200000 | 01222fttt | tttt = ttt.t degrees f = 0 Celsius f = 2 Fahrenheit |
| Print Frequency | Read | 0123200000 | 0123200mmm | mmm = 0-255 min |
| | Write | 0133200mmm | Same | |
| Status | Read | 0125000000 | 01250vltep | v = VK 7000 control l = low level t = over temp. e = ext. sensor p = printer on 1/0 = Yes/No |
| Print Instantly | Write | 0135100000 | Same | |
| Reset | Write | 0135000000 | Same | |

RS232 Connections

Both a standard DB-9 subminiature connector and RJ-11 modular connectors are provided for digital communications with an external controller such as a PC. The pertinent pin connections are detailed in Figure 7, “DB-9 Connector for RS232,” below.

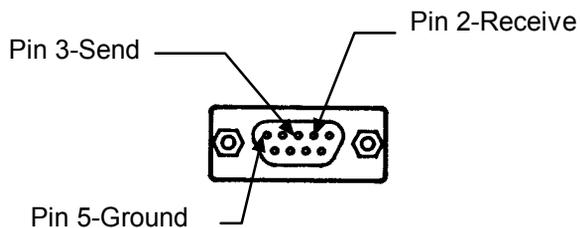
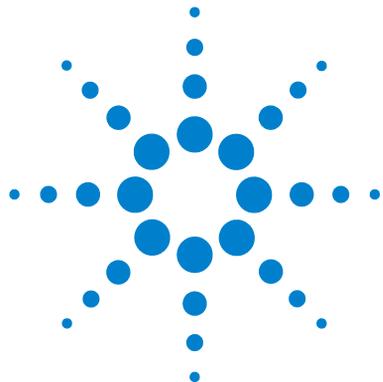


Figure 7 DB-9 Connector for RS232

Control software, which runs in Microsoft Windows®, is available from the Dissolution Systems Service Department. Please contact the Dissolution Systems Service Department or your local Agilent representative for more information.

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



5 Maintenance

Periodic Maintenance 42



Periodic Maintenance

WARNING

The instrument 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.

Periodic maintenance intervals may vary depending on frequency of instrument usage.

Monthly

- Clean and dry all surfaces.
- Apply an algaecide, bactericide, or any other additive that will not corrode plastic or stainless steel to the water bath or system.
- If the water is high in calcium, a 10% solution of white vinegar and water can be used to remove deposits. *Do not use this solution more than once a month.*
- Flush external tubing and check clamps for corrosion.
- Clean the external temperature probe with alcohol to remove deposits. Deposits can reduce the measurement accuracy.
- If a four-pin communication cable is used, inspect the cable for damage or corrosion. Inspect the power entry module for corrosion and clean it if necessary. Replace the fuse as necessary.

Fuse Replacement

CAUTION

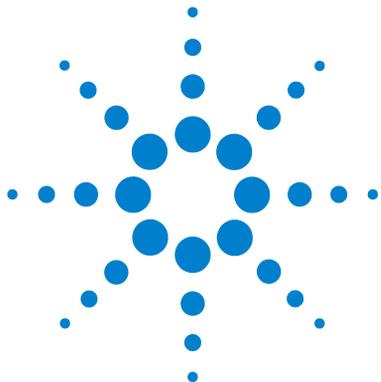
Panels or covers that are retained by fasteners which require the use of a tool for removal may be opened only by Agilent-trained, Agilent-qualified, or Agilent-authorized service engineers.

The line fuse is located in a small cassette built into the rear panel AC power receptacle. To remove the line fuse, press the release tab down and slide the entire cassette out of the receptacle. Snap the fuse out of the cassette, replace the fuse, and reinsert the cassette pushing until the release tab snaps back into place.

WARNING

Never replace a fuse with a higher rated fuse. Use *only* a 10 amp fuse for 115 V power or *only* a 5 amp fuse for 230 V.

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



6 Service and Warranty

Service and Warranty Information 46

The warranty is provided by Agilent Technologies, Inc. or one of its authorized representatives.



Service and Warranty Information

Agilent dissolution products carry a one-year warranty on parts and labor. The Dissolution Systems Service Department (or one of its representatives) will, at its option, either repair or replace any mechanical and electrical components in your instrument which prove to be defective. During the first year of warranty coverage, there is no charge for the labor to repair your unit. The Dissolution Systems Service Department (or one of its representatives) will determine the best site to repair the unit, either onsite or returned to Agilent Technologies, Inc. Any onsite warranty services are provided only at the initial installation point. Installation and onsite warranty services are available only in Dissolution Systems service travel areas.

Exclusions and Limitations

Excluded from this warranty are expendable or consumable items such as, but not limited to, paddles, baskets, vessels, and acrylic water baths. Also excluded are defects from improper or inadequate maintenance by the customer, user-induced chemical action or contamination, unauthorized modification or misuse, and improper site preparation and maintenance.

Operation of software is not warranted to be uninterrupted or error-free.

Obtaining Warranty Service

To obtain warranty service in the United States, contact the Dissolution Systems Service Department at 800.229.1108 to obtain authorization to return units for repair. At the option of the customer, onsite warranty service is available, but travel charges may be incurred. The customer should prepay all shipping charges for products returned to the Dissolution Systems Service Department (unless otherwise authorized), and Agilent Technologies, Inc. will pay all charges for return to the customer.

Warranty Limitations

Agilent Technologies, Inc. makes no other warranty, either express or implied, with respect to this product. Specifically disclaimed are any implied warranties of merchantability and fitness for a particular use. In no event will Agilent Technologies, Inc. be liable for any indirect, incidental, or consequential damages arising from the use of this product. This warranty gives you specific legal rights which may vary from state to state or province to province, so you may have other rights and some of these exclusions may not apply to you.

Exclusive Remedies

The remedies provided herein are the customer's sole and exclusive remedies. In no event shall Agilent Technologies, Inc. or its representatives be liable for any direct, indirect, special, incidental, or consequential damages, whether based on contract, tort, or any other legal theory. Some states or provinces do not allow the exclusion or limitation of incidental or consequential damages, so the above limitation or exclusion may not apply to you.

6 Service and Warranty

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

Index

Numerics

750D Keypad, 34
750D Rear Panel, 21

A

Apparatus 1 / 2 Control, 32
Apparatus 3 / Apparatus 7
Control, 33

C

Conventions, 17

D

DB-9 Connector for RS232, 39
Dual Water Bath Tubing
Connections, 25

E

Electrical Connections, 21
Exclusions, 46
Exclusive Remedies, 47

H

Hidden Key Functions, 27

I

Initial Power Up, 30
Interfacing to the dissolution
apparatus, 22

L

Limitations, 46
Liquid Connections, 23, 25

Local Control, 34

M

monthly maintenance, 42

O

Obtaining Warranty Service, 46
operating parameter values, 15
Over Temp, 35

P

Periodic Maintenance, 42
Priming the System, 31

R

Report Center Printer, 36
RJ-11 modular connectors, 39
RS232 Command Set, 38
RS232 Connections, 39

S

Service, 46
Single Water Bath Connections, 23

U

Unpacking, 20

W

Warranty, 46
Warranty Limitations, 47

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