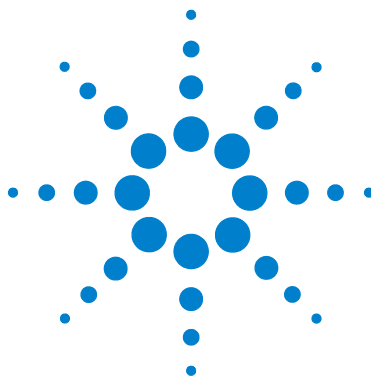




Rotating Bottle Apparatus

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

70-9062

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
1 Safety	7
Electrical Hazards	7
Warning	9
Caution	9
Note	9
Information Symbols	10
General	11
WEEE Directive	12
2 Introduction	13
Conventions Used in this Manual	15
3 Setting Up the Rotating Bottle Apparatus	17
Unpacking Your Equipment	18
Setting up the Water Bath and 750	19
Connecting the Temperature Probe	20
Initial Power Up	20
Power Cord Connections	20
Powering Up the 750 Heater / Circulator	21
Powering Up the Rotating Bottle Apparatus	21
4 Operating the Rotating Bottle Apparatus	23
Operating Instructions	24

Contents

5	Troubleshooting and Maintenance	27
	Periodic Maintenance	28
	Daily Maintenance	28
	Weekly Maintenance	28
	Monthly Maintenance	29
	Annual Maintenance	29
	Water Bath / Acrylic Care	30
	Fuse Replacement	31
	Troubleshooting	32
6	Service and Warranty	33
	Service and Warranty Information	34
	Exclusions and Limitations	34
	Obtaining Warranty Service	34
	Warranty Limitations	35
	Exclusive Remedies	35
	Index	37

Figures

Figure 1. Rotating Bottle Apparatus [14](#)

Figures

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



1 Safety

Electrical Hazards	7
WEEE Directive	12

The Rotating Bottle Apparatus 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 Rotating Bottle Apparatus involves the use of solid dosage forms and 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 Rotating Bottle 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 8020 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 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 Rotating Bottle Apparatus 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 Rotating Bottle Apparatus 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 wheelie bin” WEEE symbol shown on page 10 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 15

The Rotating Bottle Apparatus is designed for release rate testing of tablets or capsules. Manual changing of media during a test simulates a dosage passing through a pH gradient.

The Rotating Bottle Apparatus includes the following items:

- One system drive with easy-to-read digital display
- One clear acrylic water bath, 19 x 12.7 x 15.3 inches, with 750 Heater / Circulator or one clear acrylic water bath, 24 x 12 x 12 inches, for use with an immersion circulator
- One stainless steel drive bar with 14 positions
- 20 standard 100 mL test bottles with caps
- Five decant caps with 40-mesh screen

WARNING

The Rotating Bottle 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



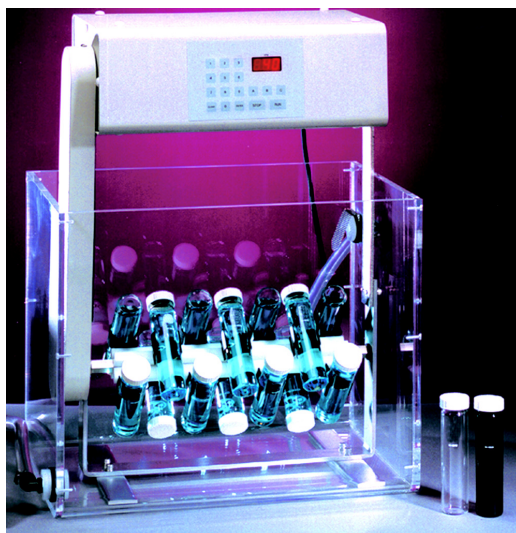


Figure 1 Rotating Bottle Apparatus

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”.

NOTE

Remember to return the warranty card supplied with this manual. 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 that arise concerning your equipment, such as upgrades, retrofits, or regulatory changes.

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



3

Setting Up the Rotating Bottle Apparatus

Unpacking Your Equipment 18

Setting up the Water Bath and 750 19

Initial Power Up 20

WARNING

Keep your hands clear of the apparatus as it rotates.



Unpacking Your Equipment

- 1 Open each carton and check the contents for damage which may have occurred during shipping. Shipping damage rarely occurs, but if it does contact both the carrier who delivered the instruments and the Dissolution Systems Service Department. Though claims for damage should be filed with the carrier, we can help you file a claim.
- 2 Carefully remove the Rotating Bottle Apparatus from its shipping carton. Hold the unit firmly to prevent it from dropping.

WARNING

Hold the apparatus only by the base. To avoid damage, do not hold the apparatus by the drive unit.

- 3 Check for items which may have come loose during shipping. Be sure to remove all parts before discarding or storing the packaging.
- 4 Place the unit on a clean, dry, level section of the bench top or table. At least four inches (10 cm) of space should be open at the rear and on both sides of the instrument.

WARNING

The electrical connection at the back of the apparatus is the primary disconnect for the instrument. The apparatus should be positioned to allow accessibility to the power cord for easy disconnection.

Setting up the Water Bath and 750

NOTE

If your equipment is not supplied with a 750, attach the heater / circulator according to the manufacturer instructions.

Complete the following steps to install the 750:

- 1 Locate the two lengths of 1/2-inch plastic tubing (one long, one short) and four stainless steel tubing clamps in the 750 tubing kit.

NOTE

These instructions assume the 750 is placed to the right of the Apparatus 7 when looking at the front of the equipment. If your 750 is placed to the left of the Apparatus 7, use the longer tubing when the instructions call for the shorter tubing and the shorter tubing when the instructions call for the longer tubing.

- 2 Slide a tubing clamp over one end of the shorter length of tubing supplied. Carefully place this end of the tubing over the connector labeled INLET on the 750 rear panel. Slide the tubing on until it meets the rear panel.
- 3 Slide the tubing clamp over the connector and tighten securely.
- 4 Place a tubing clamp over the opposite end of the same piece of tubing. Attach the tubing to the open end of the Y connector coming from the barbed-angle tubing adapter on the right side of the water bath.
- 5 Slide the tubing clamp over the connector and tighten securely.
- 6 Connect the water bath inlet located on the rear of the water bath to the 750 rear panel connector labeled OUTLET in the same fashion using the longer piece of tubing.
- 7 Connect the four-pin DIN cable between the jack labeled INPUT SIGNAL on the 750 rear panel and the jack labeled HEATER CIRCL. on the Apparatus 7 right side panel.

- 8 Ensure the 750 rear panel power switch is in the OFF position.

NOTE

If possible, preheat the water to speed achievement of the desired water bath temperature. Use purified water whenever possible to minimize scale and mineral buildup. Algaecides may be used to inhibit algae and bacteria growth. Check the label to ensure the formulation is compatible with the plastic materials used in the water bath construction.

Connecting the Temperature Probe

- 1 Plug the supplied water bath temperature probe into the jack labeled BATH TEMP on the Apparatus 7 right side panel.
- 2 Place the temperature probe through the small hole in the vessel table on the rear left side so that approximately 3/4 of the probe is below the surface of the table.

Initial Power Up

Power Cord Connections

- 1 Ensure the Rotating Bottle Apparatus is turned off.

WARNING

Before plugging the Rotating Bottle Apparatus into any power outlet, ensure the instrument is properly configured for the voltage provided. Check the serial number tag on the rear panel of the instrument to confirm the voltage requirement.

The electrical connection at the back of the apparatus is the primary disconnect for the instrument.

- 2 Connect power cords to the receptacles at the rear of the Rotating Bottle Apparatus and the 750.
- 3 Plug the cords into electrical outlets of the appropriate voltage.

Powering Up the 750 Heater / Circulator

Turn on the 750. Flow should begin immediately and the unit should operate quietly with a low “hum”. Bubbles may appear at the water bath inlet as air is expelled from the system. After a few minutes flow into the water bath should be smooth and steady. If not, stop the system and prime it as described in the *750 Heater / Circulator Operator’s Manual*.

NOTE

Make a final check for leaks. Allow the heater / circulator to run for five minutes, then check the water bath, tubing, connectors, and the bench area immediately surrounding the apparatus. If any water leaks are seen, turn off the power immediately and remove the power cord. Locate and fix the source of the leak before proceeding.

Powering Up the Rotating Bottle Apparatus

Turn on the Rotating Bottle Apparatus. The red LED on the front panel illuminates. If it does not, ensure the unit is securely plugged in and there is power at the outlet. If the unit still fails to respond, see [“Troubleshooting”](#) on page 32 or call the Dissolution Systems Service Department.

3 Setting Up the Rotating Bottle Apparatus

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



4 Operating the Rotating Bottle Apparatus

Operating Instructions 24

WARNING

Keep fingers and loose clothing away from the lower end of the drive belt assembly.



Operating Instructions

The procedure as described in NF XIV is not an official method, but is suggested as a suitable test that could be helpful in assuring product uniformity of timed-release tablets and capsules. See General Tests, “Timed-Release Tablets and Capsules—In Vitro Test Procedure,” pages 985-986, NF XIV, July 1, 1975.

To operate the Rotating Bottle Apparatus, complete the following steps:

- 1 Add tablets or capsules and extracting fluid to the standard test bottles.
- 2 Secure the standard caps on the bottles.
- 3 Snap the bottles into the clips.

NOTE

For even speed control, the bottles should be distributed in parallel pairs on opposite sides of the shaft.

- 4 Center each bottle within the clip.
- 5 Carefully place the apparatus in the water bath.
- 6 Turn on the Rotating Bottle Apparatus.
- 7 Press **C** on the keypad to select the speed and enter the desired rotation speed (between 10 RPM and 60 RPM).

Optionally, press **B** on the keypad to toggle between 30 RPM and 40 RPM.

NOTE

Key A on the keypad is inactive.

- 8 Press **RUN** to begin rotation of the bottles.
- 9 At the appropriate sampling time, press **STOP**. The apparatus pauses and the bottles stop rotating.
- 10 Remove the Rotating Bottle Apparatus from the water bath.
- 11 Remove the bottles from the clips.
- 12 Replace the standard caps with 40-mesh decant screen caps.

- 13 Decant, retaining the residue.
- 14 Add extracting fluid to the residue.
- 15 Replace the 40-mesh decant screen caps with the standard caps.
Ensure the caps are secured tightly.
- 16 Snap the bottles into the clips.

NOTE

For even speed control, the bottles should be distributed in parallel pairs on opposite sides of the shaft.

- 17 Center each bottle within the clip.
- 18 Carefully place the Rotating Bottle Apparatus in the water bath.
- 19 Press **RUN** to continue the test.
- 20 When the test is complete, press **STOP**.

4 Operating the Rotating Bottle Apparatus

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



5 Troubleshooting and Maintenance

Periodic Maintenance 28

Troubleshooting 32



Periodic Maintenance

WARNING

The Rotating Bottle 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.

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

Daily Maintenance

- Immediately wipe all spills or errant materials from the exterior of the instrument or any exposed part using a clean cloth and plain water.
- Carefully wash the bottles and caps after each use.

Weekly Maintenance

See “[Water Bath / Acrylic Care](#)” on page 30 as applicable for additional information on proper maintenance of your equipment.

- Inspect the water bath and heater / circulator tubing for algae or other materials. If algae is present, change the bath water and add an algicide.
- If you use a water bath algicide or clear bath product, ensure it is compatible with PETG and acrylic.

Monthly Maintenance

See “Water Bath / Acrylic Care” on page 30 as applicable for additional information on proper maintenance of your equipment.

- Clean the outside of the apparatus with a damp cloth.
- Drain the water from the water bath and clean the bath thoroughly. Refill the water bath and add an algicide.
- It is recommended that the water bath temperature probe jack is checked for surface corrosion and wiped clean with a soft cloth or nonabrasive pad every one to three months.

NOTE

Depending on the frequency of use, it may be necessary to complete this maintenance procedure more often.

Annual Maintenance

- Inspect all electrical connections for corrosion and damage.
- Clean the outside of the apparatus with a damp cloth.
- Check the belt tension—total deflection should not exceed 1/2 inch (1.3 cm).

CAUTION

Do not use cleaning compounds containing ammonia or abrasive cleaners on your water bath.

Water Bath / Acrylic Care

The water bath supplied with the Rotating Bottle Apparatus should be maintenance free except for an occasional cleaning. If you use a water bath algaecide or clear bath product, ensure it is compatible with PETG and acrylic. The flow paths in the immersion circulator are primarily stainless steel and should tolerate most clear bath formulations. Check with the product manufacturer to be sure the product is safe for your water bath.

- All of our water baths are fabricated entirely of commercial grade acrylic. When using them 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 PETG and acrylic before use. If in doubt, call the service department for advice before proceeding.
- Do not use ammonia, window-cleaning sprays, kitchen scouring compounds, or solvents such as acetone, gasoline, benzene, alcohol, carbon tetrachloride, or lacquer thinner. These can scratch the material's surface and / or weaken it causing small surface cracks called "crazing".

- Our recommendations include but are not limited to the following:
 - Hot water: < 150 °F
 - Vinegar (5% Glacial Acetic Acid)
 - Ethyl alcohol: maximum 10%
 - Isopropyl alcohol: maximum 25%

Fuse Replacement

WARNING

The Rotating Bottle 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.

Follow these steps to replace the fuse:

- 1 Before checking or attempting to replace a fuse, remove the power cord from the unit.
- 2 Press the release tab on the fuse holder located in the power entry module with the blade of a small screwdriver or equivalent.
- 3 Upon release, slide the holder out of the AC power entry module.
- 4 Remove the old fuses and insert the new fuses.
- 5 Slide the holder back into the power entry module and push until it locks into place.
- 6 Replace the power cord.

WARNING

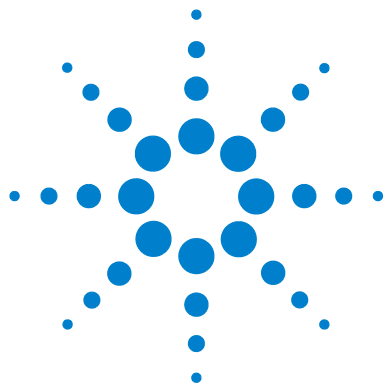
Never replace a fuse with one of a higher amperage rating. Doing so may compromise the safety margin and could result in damage to the instrument or personal injury.

Troubleshooting

The Dissolution Systems Service Department can assist you if you experience problems or have questions concerning your Rotating Bottle Apparatus. Many problems can be traced to simple sources and are easily solved.

Following is a troubleshooting guide which may help you. The Dissolution Systems Service Department can be reached at 800.229.1108 (inside the US) or 919.677.1108 (outside the US). Optionally, you can send a fax to 919.677.1138. You can also e-mail the Dissolution Systems Service Department at dissolution.service@agilent.com.

Symptom	Probable Cause	Possible Solution
There is no power to the apparatus.	The fuses are blown (open).	Replace the fuses in the power entry module. See “Fuse Replacement” on page 31.
	There is no power at the electrical outlet.	Verify that the outlet has power going to it.
	The power cord is damaged or is not plugged in.	Check the power cord for abrasion and proper connection. Replace if necessary.
	The power entry module is faulty.	Check the module to ensure proper voltage delivery. Contact the Dissolution Systems Service Department for assistance.



6 Service and Warranty

Service and Warranty Information 34

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.

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

Index

A

Acrylic care, [30](#)
Annual Maintenance, [29](#)

C

Conventions, [15](#)

D

Daily Maintenance, [28](#)

E

equipment, unpacking, [18](#)
Exclusions, [34](#)
Exclusive Remedies, [35](#)

F

Fuse Replacement, [31](#)

H

Heater / Circulator, powering up, [21](#)

L

Limitations, [34](#)

M

Monthly Maintenance, [29](#)

N

No power, [32](#)

O

Obtaining Warranty Service, [34](#)

P

Periodic Maintenance, [28](#)
Photograph, [14](#)
Powering Up, Heater / Circulator, [21](#)

S

Service, [34](#)

T

Temperature Probe Connection, [20](#)
Troubleshooting, [32](#)

U

unpacking your equipment, [17](#), [18](#)

W

Warranty, [34](#)
Warranty Limitations, [35](#)
Water bath Care, [30](#)
Weekly Maintenance, [28](#)

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