

100 Automated Disintegration Apparatus

Operator's Manual



Notices

Manual Part Number

SD-29001863 Rev. E
Edition: 03/2022

Copyright

© Agilent Technologies, Inc. 2020, 2022

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

This manual covers the following modules:

- Agilent 100 Automated Disintegration Apparatus (G7962A)

1 Safety 6

Electrical Hazards 7

Warning 8

Caution 8

Note 8

Information Symbols 9

2 Introduction 10

Introduction 11

Conventions Used in this Manual 12

Serial Number Format 12

3 Setting Up the 100 13

Unpacking Your 100 14

Environmental Requirements for Installation 15

Setting up the Water Bath 16

Initial Power Up 17

Electrical Connections 17

Content

Installing the Temperature Probe	19
Water Bath Status LEDs	20
Setting the Date / Time Clock	21
Mounting Baskets and Filling Beakers	22
Draining the Water Bath	22
Hidden Key Functions	23
4 Operating the 100	24
Setting Water Bath Temperature	25
Setting Print Frequency for the Report Center Printer	25
Adjusting Basket Height	26
Placing Dosage Units in the Baskets	27
Selecting Operating Mode	27
Single-basket Operation	27
Multi-basket Operation	28
Three-Basket Operation	31
Printer Operation During Manual or Program Operation (Optional)	31
5 Troubleshooting and Maintenance	32
Preventive Maintenance	33
Weekly Maintenance	33
Monthly Maintenance	33
Every Six Months Maintenance	34
Yearly Maintenance	34
Disintegration Basket Care	35
Water Bath / Acrylic Care	36
Visual Checks	37
Water Bath Temperature Probe Accuracy Test	38

Content

Report Center Impact Printer 40

Installing the Cartridge Ribbon 40

Replacing the Paper Roll 41

Toggling Your Printer Online 42

Printer Self Test 43

Printer Configuration 44

Troubleshooting 46

Obtaining Warranty and Other Services 47

1

Safety

Electrical Hazards 7

The Agilent 100 (G7962A) 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.

The Agilent 100 (G7962A) is operated in conjunction with equipment that uses 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 Agilent 100 (G7962A) 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 100 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



Switches main power on



Switches main power off



Indicates single-phase alternating current



The apparatus is marked with this symbol when the user shall refer to the instruction manual in order to protect risk of harm to the operator and to protect the apparatus against damage.



Confirms that a manufactured product complies with all applicable European Community directives. The European Declaration of Conformity is available at:

<http://regulations.corporate.agilent.com/DoC/search.htm>



Confirms that a manufactured product complies with all applicable UK Regulation.

The UK Declaration of Conformity is available at:

<http://regulations.corporate.agilent.com/DoC/search.htm>



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



This product complies with the European WEEE Directive marking requirements. The affixed label indicates that you must not discard this electrical/electronic product in domestic household waste.



NOTE

Do not dispose of in domestic household waste!

To return unwanted products, contact your local Agilent office, or see <http://www.agilent.com> for more information.



ISM1-A

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

ISM classification according to CISPR11/EN55011: ISM Group 1 Class A

CAN ICES-001(A)

NMB-001(A)

This ISM device complies with Canadian ICES-001. Cet appareil ISM est conforme à la norme NMB-001 du Canada



2

Introduction

Introduction	11
Conventions Used in this Manual	12
Serial Number Format	12

Introduction

The 100 (G7962A) is designed to provide versatility and reliability, while meeting all current USP testing requirements. Its three basket positions can be programmed to dip together or independently, allowing up to three different tests to be run at the same time. In addition to the standard six-tube USP-approved basket rack, a variety of optional basket racks are available for large and bolus tablets. The 100 is used for disintegration testing in accordance with current USP requirements. The optional built-in Report Center Printer provides hard-copy documentation of the instrument operation for the highest degree of confidence.

The apparatus is fully programmable by simply entering the time duration of the test. Baskets automatically lift from the beaker at the end of the test in the pass / fail mode. The 100 also has individual digital time displays for each basket and the water bath temperature is displayed continuously on the front panel.

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. See “Obtaining Warranty and Other Services” on page 47..

Conventions Used in this Manual

- Items you are asked to press are in bold. For example, “press **H** on the keypad”.

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 100

Unpacking Your 100	14
Setting up the Water Bath	16
Initial Power Up	17
Installing the Temperature Probe	19
Water Bath Status LEDs	20
Setting the Date / Time Clock	21
Mounting Baskets and Filling Beakers	22
Draining the Water Bath	22
Hidden Key Functions	23

Unpacking Your 100

- 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 100 from its shipping carton. Hold the unit firmly to prevent it from dropping.

WARNING

Hold the unit only by the base. Do not hold by the reciprocating shafts to avoid damage.

- 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 of space should be open for the rear panel power connections and for access to the rear panel power switch.

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 cords for easy disconnection.

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
- Overvoltage Category: II
- Altitude: 2000m
- Temperature: 5 °C to 40 °C
- Power: 115 V, 50/60 Hz, 10 A or 230 V 50/60 Hz, 5 A

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

Setting up the Water Bath

- 1 Ensure the 100 is not plugged in.
- 2 Place the water bath on the 100 with the lower quick-connect fitting positioned on the left side of the apparatus. Use the locator holes on the water bath and the locator pins on the 100 to ensure proper placement.
- 3 Locate the two tubes with quick-connect fittings and insert them into the appropriate quick-connect fittings on the 100 water bath (see Figure 1, “Water Bath Connections,” below).



Figure 1. Water Bath Connections

- 4 Ensure the cover is in place on the water bath with the small hole toward the front. Ensure the disintegration beakers are not in place in the cover.
- 5 Fill the water bath through any one of the three beaker holes in the cover.

NOTE

Initially fill the water bath no higher than 2.5 inches (6.4 cm) from the top, to allow for displacement by the beakers when they are inserted.

- 6 Check the water bath and all connecting tubing for leaks. When satisfied that no leaks are present, proceed to **“Installing the Temperature Probe”** on page 19.

Initial Power Up

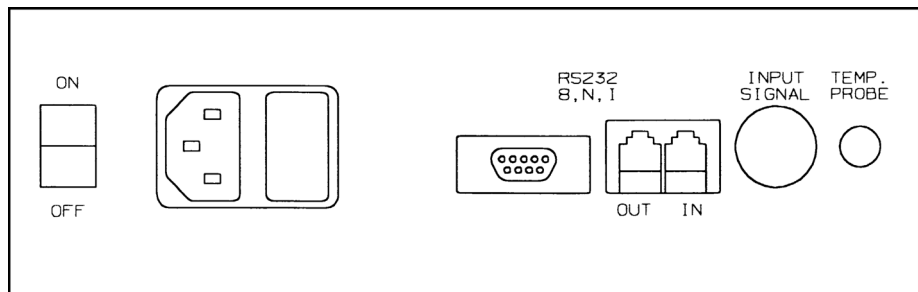
Electrical Connections

WARNING

Before plugging the 100 into any power outlet, ensure the instrument is configured for the voltage provided. Check the serial number tag on the rear panel to confirm the voltage requirement.

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

See **Figure 2**, “100 Rear Panel,” on page 17 to complete the steps that follow.



Port	Connector	Max In / Out Voltage	Function
RS 232	RS 232	12 Vdc	Communication port to PC for remote control
OUT	RJ11	n/a	Not supported
IN	RJ11	n/a	Not supported
Input Signal	4-pin DIN	n/a	Not supported
Temp Probe	6.5mm coax	5 Vdc	Probe temperature from water tank

Figure 2. 100 Rear Panel

- 1 Ensure the power is off.
- 2 Connect the power cord between the AC power connector and an AC power receptacle of the appropriate voltage.

- 3 Turn on the 100. The three position screens and the temperature screen on the front panel illuminates immediately. You will hear the circulator working as it begins to pump the bath water through the system. Bubbles may appear at the water bath inlet as air is expelled from the system.

NOTE

Check for leaks. Allow the circulator to run for five minutes, then check the 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.

- 4 Place the disintegration beaker(s) through the holes in the top cover of the water bath.
- 5 As the water circulates, there may be a slight drop in the water bath level as the internal tubing fills. Wait until the beakers are in place before adding additional water as required to bring the level back to within no more than 2.5 inches of the top of the water bath. Replace the cover, if removed.

Installing the Temperature Probe

The 100 is shipped with a factory-installed temperature probe inside the heater /circulator tank and an external probe to be placed directly in the water bath. Both probes are necessary and the temperature readings must remain within 5 °C of each other or the 100 will not operate. If either probe is not functioning or not installed, the alarm sounds and the keypad is inactive.

NOTE

No installation or positioning is necessary for the internal temperature probe.

To install the external temperature probe, complete the following steps:

- 1 Plug the temperature probe into the jack labelled TEMP. PROBE on the rear panel.
- 2 Insert the metal tip of the probe into the small hole in the water bath cover.
- 3 Ensure the small hole and temperature probe are toward the front of the water bath.
- 4 Press **INT/EXT** to view the temperature readings for each probe. When the external probe reading displays, the Ext Temp Probe LED illuminates.

Water Bath Status LEDs

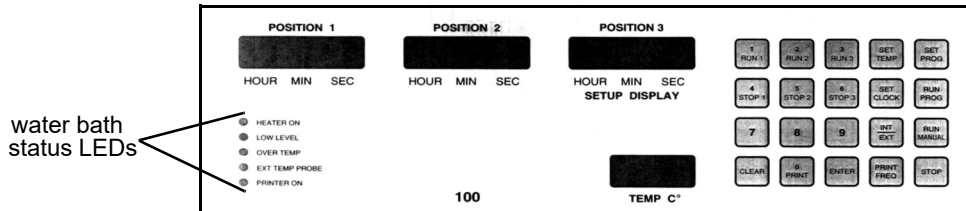


Figure 3. 100 Keypad

The Heater On LED flashes to indicate heater activity. When the heater is on continuously, as will occur when the water bath is cold, the LED remains illuminated. As the target temperature is approached, the LED flashes indicating the heater is cycling on and off.

The Over Temp LED illuminates

- if the temperature of the water of the water bath exceeds 65 °C. Check for obstructions and low water if this condition occurs. The over temp condition clears automatically when the temperature cools below 55 °C. Press **CLEAR** to reset if the temperature is between 55 and 65 °C.
- if the external temperature probe is not in the water bath or is not connected.
- if the temperature probe readings are more than 5 °C higher than the set temperature.

The Ext Temp Probe LED illuminates when the external temperature probe reading displays in the temperature screen. Otherwise the internal probe reading displays.

NOTE

If the external temperature probe is not connected, the temperature display reads 4.4. This is normal and changes as soon as the probe is installed.

The Printer On LED illuminates when the Report Center Printer has been enabled.

Setting the Date / Time Clock

If you are using the built-in Report Center Printer to document the test conditions, you must set the correct date and time. This information displays on all printouts.

To set the date and time, complete the following steps:

- 1 Press **SET CLOCK**. The Position 3 screen flashes the date.
- 2 If the date is correct, press **ENTER**. Otherwise, enter the correct date in mm/dd/yy format and press **ENTER**. The time flashes.
- 3 If the time is correct, press **ENTER**. Otherwise, enter the correct time in 24-hour format and press **ENTER**.

Mounting Baskets and Filling Beakers

- 1 Ensure the water bath cover is in place on the water bath.
- 2 Place one, two, or three 1000 mL USP disintegration beakers through the water bath cover holes so they are suspended in the water bath.
- 3 Hold the beaker(s) down and pour 900 mL of the required disintegration medium in each beaker. When full, the beakers sit securely in the cover and do not float.
- 4 Place one, two, or three basket racks onto the horizontal basket supports.
- 5 Verify the water level in the water bath is above the media level of all beakers. If not, either add additional water or drain excess water until the proper working level is reached.

Draining the Water Bath

WARNING

Disconnect the 100 from the AC power receptacle before draining the tank.

- 1 Disengage the quick-connect fitting on the left side of the tank and attach the supplied drain hose.
- 2 Open the drain valve and drain the water bath into an appropriate receptacle or sink.
- 3 Use a clean paper towel or damp cloth to wipe out the water bath and remove any remaining water from the bottom of the tank.

Hidden Key Functions

The following seldom-used operations can be performed by pressing and holding **CLEAR** and then pressing the **number** on the keypad and releasing both keys at the same time. Refer to **Figure 3**, “100 Keypad,” on page 20.

Table 1 Hidden Key Functions

Key Sequence	Function
CLEAR > 0	Use this key sequence to set the communication port identification numbers. Enter the identification number and press ENTER .
CLEAR > 2	Use this key sequence to enable or disable alarm function.
CLEAR > 5	Use this key sequence to toggle between 1 and 2 decimal places for the temperature display. Two decimal places is used to calibrate the temperature.
CLEAR > 8	Use this key sequence to select the baud rate. This key combination toggles among 1200, 2400, 4800, and 9600. When the correct baud rate displays, press ENTER .



4

Operating the 100

Setting Water Bath Temperature 25

Setting Print Frequency for the Report Center Printer 25

Adjusting Basket Height 26

Placing Dosage Units in the Baskets 27

Selecting Operating Mode 27

Printer Operation During Manual or Program Operation (Optional) 31

Setting Water Bath Temperature

To set the water bath temperature, press **SET TEMP**. Enter the desired water bath temperature and press **ENTER**.

NOTE

It may be necessary to set the water bath temperature up to 0.5 °C higher than the desired temperature of the medium in the beakers to compensate for heat loss through the walls of the beakers.

Setting Print Frequency for the Report Center Printer

If your 100 is equipped with a Report Center Printer and you want an automatic periodic printout of test conditions, enable the built-in Report Center Printer and indicate the desired automatic print interval.

To enable the printer, press **PRINT FREQ**. Enter the desired auto-print interval in mmm format up to 255 minutes and press **ENTER**.

To disable the printer, enter an auto-print frequency of **0** (zero) and press **ENTER**.

Adjusting Basket Height

- 1 Lower the basket to the lowest position and measure the distance from the bottom of the beaker to the bottom of the basket. This measurement should be no less than 2.5 cm.
- 2 If necessary, adjust the basket height by loosening the screw on the basket hanger with an Allen wrench. Extend or retract the two parts and tighten the screw to achieve the proper distance of not less than 2.5 cm from the bottom of the beaker.

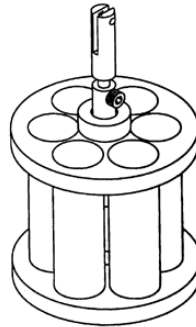


Figure 4. Basket Adjustment

Placing Dosage Units in the Baskets

Place one dosage unit in each of the six tubes of the basket. USP fluted disks may be added where specified in the individual monograph.

Selecting Operating Mode

You can select single-basket, multi-basket, or three-basket operation.

NOTE

The Position 3 screen is the display screen for all programming regardless of which position or parameter you are programming.

Single-basket Operation

In single-basket mode, the operation of each basket is set independent of the others. To start a single basket:

- 1 Press **RUN #** (where # is the basket number you want to run). The current test time for that position, if any, flashes. For example, the following display indicates that the basket in the selected position is currently programmed to dip for 10 minutes:

POSITION 3

001000

- 2 To accept the current test length, press **ENTER**. The basket lowers and begins dipping, lifting automatically at the end of the programmed time period.
To enter a different dip time value, enter the new value in hhmmss format and press **ENTER** to start the dip cycle.
- 3 To stop the dip cycle of a basket before the end of the programmed time, press **STOP #** (where # is the basket number you want to stop). The basket stops dipping and lifts from the beaker immediately.

NOTE

Press **STOP** to immediately stop *all* baskets.

Multi-basket Operation

There are two benefits of multi-basket mode:

- You can program any combination of baskets to start and stop at exactly the same time. This feature allows the 100 to function as a conventional two- or three-basket apparatus.
- You can program the 100 to drop the baskets at staggered times and to remove them at the same time. This feature is useful when running different products with different test lengths simultaneously without returning to the apparatus at different times to view the end of each test.

For example, you can position baskets 1 and 3 to start and finish their dip cycles at exactly the same time while Position 2 remains idle, or baskets 1 and 2 start and finish while Position 3 remains idle. Alternatively, you can have basket 3 start automatically 5 minutes after basket 1, with a test time 5 minutes less than that for basket 1. This causes both baskets to finish at the same time while Position 2 remains idle.

The screen under Position 3 acts as the display screen for programming the 100. Create a programming sequence as follows:

- 1 Press **SET PROG**. In the display, the numeral 1 flashes, alternating with the previously programmed test length for that position in hhmmss format. For example, you may see the following alternating displays:

POSITION 3

1

POSITION 3

003000

The example on page 28 indicates the previously programmed test length for Position 1 was 30 minutes.

- To accept the current test length, press **ENTER**.
 - To change it, enter the new value in hhmmss format and press **ENTER**.
 - To disable Position 1, enter **000000** for the test length and press **ENTER**.
- 2** In the display, the numeral 2 flashes, alternating with the previously programmed delay time for Position 2 in hhmmss format. The delay time is the amount of time between the start of the program and the start of dipping for Position 2. For example, you may see the following alternating displays:

POSITION 3

2

POSITION 3

000500

The above example indicates the previously programmed delay time for Position 2 was 5 minutes.

- To accept the current delay time, press **ENTER**.
- To change it, enter the new value in hhmmss and press **ENTER**.
- If no delay is desired, enter **000000** and press **ENTER**.

NOTE

There is no delay time for Position 1.

- 3** In the display, the numeral 2 flashes again, this time alternating with the previously programmed test length for Position 2. For example, you may see the following alternating displays:

POSITION 3

2

POSITION 3

004000

The example on page 29 indicates the previously programmed test length for Position 2 was 40 minutes.

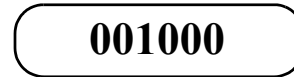
- To accept the current test length, press **ENTER**.
- To change it, enter the new value in hhmmss and press **ENTER**.

- To disable Position 2, enter **000000** and press **ENTER**.
- 4 In the display, the numeral 3 flashes, alternating with the previously programmed delay time for Position 3 in hhmmss format. The delay time is the amount of time between the start of the program and the start of dipping for Position 3. For example, you may see the following alternating displays:

POSITION 3



POSITION 3



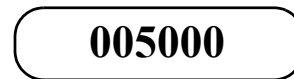
The above example indicates the previously programmed delay time for Position 3 was 10 minutes.

- To accept the current delay time, press **ENTER**.
 - To change it, enter the new value in hhmmss and press **ENTER**.
 - If no delay is desired, enter **000000** and press **ENTER**.
- 5 In the display, the numeral 3 flashes again, this time alternating with the previously programmed test length for Position 3. For example, you may see the following alternating displays:

POSITION 3



POSITION 3



The above example indicates the previously programmed test length for Position 3 was 50 minutes.

- To accept the current test length, press **ENTER**.
 - To change it, enter the new value in hhmmss and press **ENTER**.
 - To disable Position 3, enter **000000** and press **ENTER**.
- 6 Press **RUN PROG**. The programmed sequence displays in the Position 3 screen. If correct, press **ENTER**. The program begins.

To stop any individual basket during multi-basket operation, press **STOP #** (where # is the basket number you want to stop). To stop all baskets at the same time, press **STOP** in the lower right corner of the keypad.

NOTE

When an individual basket is stopped and the other baskets are dipping, the individual basket can only be restarted in the independent single-basket mode. It cannot be restarted as part of the running program.

Three-Basket Operation

To run all three baskets together, starting and stopping at the same times, complete the following steps:

- 1 Press **RUN MANUAL**. The previously set test length (if any) for three-basket operation flashes in the Position 3 screen.
- 2 Press **ENTER** to accept the displayed values or enter a new test length in hhmss format and press **ENTER**. All three baskets lower and begin dipping. The baskets stop simultaneously, lifting automatically at the end of the programmed time period.

NOTE

Press **STOP** to stop all three baskets at any time during the test.

Printer Operation During Manual or Program Operation (Optional)

If your 100 is equipped with the optional Report Center Printer, you can generate a printout of the instrument status at any time by pressing **PRINT**. This does not affect automatic printing as set by the PRINT FREQ function.



5

Troubleshooting and Maintenance

Preventive Maintenance 33

Water Bath Temperature Probe Accuracy Test 38

Report Center Impact Printer 40

Troubleshooting 46

Preventive Maintenance

WARNING

The 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.

Preventive maintenance intervals may vary depending on frequency of instrument usage. To place an order for warranty or other services, see **“Obtaining Warranty and Other Services”** on page 47.

Weekly Maintenance

See **“Disintegration Basket Care”** on page 35 and **“Water Bath / Acrylic Care”** on page 36 as applicable for additional information on proper maintenance of your equipment.

- Ensure media is removed and the system is wiped with a clean, damp cloth. Inspect the water bath and circulator tubing for algae, other material and cracks / damage. If algae is present, change the bath water and add algaecide.
- If any stainless steel parts show signs of surface discoloration, lightly wipe the surface with a soft cloth or nonabrasive pad to remove it.

Monthly Maintenance

See **“Water Bath / Acrylic Care”** on page 36 as applicable for additional information on proper maintenance of your equipment.

- Ensure media is removed and the system is wiped with a clean, damp cloth.
- Lightly spray Lube1/ WD-40™ onto the shafts and run the machine for a few minutes.
- Change the bath water and add algaecide.

Every Six Months Maintenance

See **“Water Bath / Acrylic Care”** on page 36 as applicable for additional information on proper maintenance of your equipment.

- Ensure media is removed and the system is wiped with a clean, damp cloth.
- Lightly spray Lube1/ WD-40™ onto the shafts and run the machine for a few minutes.
- Change the bath water and add algaecide.
- Inspect the tubing for algae, foreign material, and cracks / damage.

Yearly Maintenance

See **“Water Bath / Acrylic Care”** on page 36 as applicable for additional information on proper maintenance of your equipment.

- Ensure media is removed and the system is wiped with a clean, damp cloth.
- Lightly spray Lube1/ WD-40™ onto the shafts and run the machine for a few minutes.
- Change the bath water and add algaecide.
- Inspect the tubing for algae, foreign material, and cracks / damage.
- Inspect the security of wires and wiring harnesses.
- Inspect the drive motor—ensure no leaking of oil has occurred.

Disintegration Basket Care

- All of our disintegration baskets are fabricated entirely of commercial grade acrylic, stainless steel, and glass. 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 acrylic before use. If in doubt, call the service department for advice before proceeding.
- Do not use 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%
- Please store disintegration baskets properly between uses. Do not simply place disintegration baskets in a drawer. They will be subject to nicks, chips, and scratches as they bump against each other. Place them back into the original shipping containers or other appropriate containers between uses. This will prevent them from coming into contact with each other or anything else in the storage area.

Water Bath / Acrylic Care

CAUTION

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

The water bath supplied with the 100 Disintegration 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 heater / 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%

Repairing Leaking Fittings

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

- 1 Turn off the heater / circulator and drain the water bath completely.
- 2 Remove the leaky bulkhead fitting.
- 3 Remove the elbow fitting from the bulkhead fitting.
- 4 Inspect the bulkhead fitting gaskets for damage and replace them as necessary.
- 5 Remove the old PTFE tape from all male fittings. Inspect the threads for damage and replace the fitting as necessary.
- 6 Apply new PTFE tape to the male fitting threads.
- 7 Reinstall and tighten the bulkhead fitting on the water bath.
- 8 Reinstall and tighten the elbow fitting to the bulkhead fitting.
- 9 Fill the water bath and turn on the heater / circulator.
- 10 Inspect the fitting for leaks. If the fitting still leaks, contact the Dissolution Systems Service Department.

Visual Checks

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. See "Obtaining Warranty and Other Services" on page 47..

Water Bath Temperature Probe Accuracy Test

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

- 1 Place the probe into a known-temperature water bath and allow several minutes for the probe to equilibrate.
- 2 The resistance is measured by attaching the leads of the voltmeter to the top and shank of the 1/4 inch phone plug.
- 3 Refer to the following table 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 ± 0.2 °C. Probes found to be outside the tolerance of ± 0.2 °C should be replaced.

Table 2 Probe Resistance Values

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

Table 2 Probe Resistance Values

Temp. Centigrade	Resistance Ohms	Temp. Centigrade	Resistance Ohms	Temp. Centigrade	Resistance Ohms
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		

Report Center Impact Printer

The following is helpful information for using your impact printer.

Installing the Cartridge Ribbon

If the printer is used infrequently, the print impression sometimes becomes weak because the ribbon dries out. If the printed material is difficult to read and you suspect this is the cause of the problem, advance to a new section of the ribbon by pressing the printer toggle switch to the **Paper feed** position. If the printing is still faint, replace the cartridge.

To install the cartridge:

- 1 Toggle the printer off line by pressing the printer toggle switch to the **OnLine / Off Line** position. When the printer is off line, the Ready LED does not illuminate.
- 2 Four small grooves are embossed on the printer cover. Gently push on these grooves to tilt the cover. When the printer cover is tilted up, you can lift it off completely.
- 3 Push down on the right side of the ribbon cartridge (marked PUSH) and remove the old cartridge.
- 4 Install the new cartridge. If there is already paper in the printer, hold the cartridge between your thumb and index finger, slide it over the paper and into the printer compartment. Ensure the paper is between the ribbon cartridge and the ink ribbon. Ensure the ink cartridge is inserted firmly to prevent weak or irregular printing. The cartridge must be properly seated and aligned for the best printing.
- 5 Turn the cartridge knob (marked by an arrow) clockwise to stretch the ribbon taut.
- 6 Replace the cover.
- 7 Toggle the printer online by pressing the printer toggle switch to the **OnLine / Off Line** position. The Ready LED illuminates.
- 8 Replace the paper if necessary.

If you get ribbon ink on the printer's plastic cover, remove it immediately. Once dried, it is difficult to remove.

Replacing the Paper Roll

- 1 Toggle the printer off line by pressing the printer toggle switch to the **OnLine / Off Line** position. When the printer is off line, the Ready LED does not illuminate.
- 2 Grasp the paper roll cover firmly by the grooves on the side and the front edge. Pull outward to remove the cover.
- 3 Press the printer toggle switch to *Paper feed* to advance the paper approximately one inch beyond the paper cutter.
- 4 Using scissors, cut the paper feeding to the printer and remove the paper roll.
- 5 Pull the remaining paper through the printer mechanism. *Pull the paper from the front (paper cutter side)*. Pulling the paper out of the back of the printer will damage the print mechanism.
- 6 Unroll several inches of paper on the new roll.
- 7 If it is jagged, cut a straight edge on the paper roll to facilitate the entry of the paper into the printer.
- 8 Slide the paper through the slot connecting the paper compartment and the printer compartment. You can slide it in approximately 1/4 inch before it stops.
- 9 While holding the paper in place, press the printer toggle switch to the *Paper feed* position and hold until approximately one inch of paper has emerged from the top of the printer.

CAUTION

Ensure the roll of paper feeds squarely. If it does not, the paper can jam and possibly damage the printer mechanism.

- 10 Release the printer toggle switch.
- 11 Turn the paper roll to take up any slack in the paper feeding to the printer.
- 12 Place the paper roll into the paper compartment.
- 13 Replace the paper roll cover. If the cover is difficult to remove or replace, the left and right edges can be trimmed or shaved with a utility knife allowing the cover to slide easier.
- 14 Toggle the printer online by pressing the printer toggle switch to the **OnLine / Off Line** position. The Ready LED illuminates.

toggling Your Printer Online

Complete these steps to toggle your printer online:

- 1 Toggle the printer online by pressing the printer toggle switch to the **OnLine / Off Line** position. When the printer is off line, the Ready LED does not illuminate.
- 2 Release the switch and it returns to the center position. The Ready LED illuminates and a READY message prints if the PRINT READY command has not been turned off. See **"Printer Configuration"** on page 44 for instructions on turning on and off the PRINT READY command. When you first turn on the instrument, it prints a READY message to assure you that the built-in microprocessor is operating properly.

When you turn off the printer, wait at least three seconds before turning it on again.

Printer Self Test

You can test the print head and ribbon only *after* inserting paper. Do not attempt to print without paper. Follow these steps to perform a printer self test:

- 1 Turn off the 100.
- 2 Press and hold the printer toggle switch in the **Paper feed** position.
- 3 Turn on the 100.
- 4 Hold the printer toggle switch until printing begins. The printer prints a list of the current configuration settings and performs a continuous print test.
- 5 Press the printer toggle switch to the **OnLine / Off Line** position to stop the printing operation.
- 6 The printer is ready to resume normal operation.

Printer Configuration

NOTE

The printer configuration is set by the factory. This procedure should be performed only if the printer displays erroneous characters. Contact the Dissolution Systems Service Department for assistance, if necessary.

- 1 Turn off the 100.
- 2 Press and hold the printer toggle switch in the **OnLine / Off Line** position while turning on the instrument. Hold the printer toggle switch in the **OnLine / Off Line** position for six seconds after the instrument is turned on, then release the switch.
- 3 The printer should print: ***** SETUP MENU ***** and **CONFIGURE... [NEXT/OK]**. If this message does not print, repeat steps 1 through 3.
- 4 The printer toggle switch is used to complete the configuration. Pressing the left side of the printer toggle switch selects **NEXT** to advance to the next menu item. Pressing the right side of the printer toggle switch selects **OK** to accept what is stated on this line of the menu item. Each time the switch is pressed, another part of the menu prints. Allow the printer to finish printing before pressing the switch again. See the table of commands on the following page.

NOTE

The printout is easier to read if the printer cover is removed.

Table 3 Printer Commands

*** SETUP MENU***		
CONFIGURE	[NEXT/OK]	Press NEXT to avoid configuration
CUSTOM	[NEXT/OK]	Press OK to enter custom mode
CUSTOM MENU		
PRINT CUSTOM SETUP	[NEXT/OK]	Press NEXT
AUTO SEQ = NO	[NEXT/OK]	Press OK
ZERO = Ø	[NEXT/OK]	Press OK
POUND SIGN = #	[NEXT/OK]	Press OK
_(UNDERScore)	[NEXT/OK]	Press OK
ONLINE/OFFLINE = YES	[NEXT/OK]	Press OK
EXT CH SET = NO	[NEXT/OK]	Press OK
PRINT READY = YES	[NEXT/OK]	Press NEXT
PRINT READY = NO	[NEXT/OK]	Press OK
READY...		

Troubleshooting

Table 4 Troubleshooting

Symptom	Probable Cause	Possible Solution
The displays do not illuminate.	The unit is not connected to an AC power source.	Connect the 100 to a power outlet.
	The fuse is blown.	Replace the fuse.
Dipping does not start immediately at an individual position.	Delay is programmed for the position in question.	Reprogram the delay.
The optional Report Center Printer does not function.	The printer is disabled.	Program a print frequency other than 000.
		Turn on the printer. The Ready LED illuminates.
The Report Center Printer functions, but nothing appears on paper.	The ribbon is exhausted.	Replace the ribbon cartridge.
The water does not circulate.	The tubing or pump is clogged.	Remove the back panel and inspect the tubing from the water bath to the pump. Remove the tubing from the pump and inspect the inlet of the pump for debris.
	The pump is faulty.	Contact the Dissolution Systems Service Department for assistance.
The OverTemp light remains on.	The water bath level is too low.	Add water to the water bath.
	Air is trapped in the tank or lines.	Attempt to prime the system by raising and lowering the front and rear of the instrument.
	The temperature probe may be in the wrong position or faulty.	Inspect the position and condition of the water bath temperature probe. Ensure the probe is plugged into the 100. Check the position of the probe within the water bath and ensure it is fully submerged. The temperature probe should only be used to measure the water bath temperature. Faulty readings occur if you use the probe in any other manner.

Table 4 Troubleshooting

Symptom	Probable Cause	Possible Solution
Operation is noisy.	Air is trapped in the tank or lines.	Attempt to prime the system by raising and lowering the front and rear of the instrument.
	The pump or motor is faulty.	Inspect the electrical connection from the pump / motor to the main PCB for damage or excessive wear. Replace the pump if necessary. Contact the Dissolution Systems Service Department for assistance.
	Plumbing connections are in the wrong configuration.	Ensure the lower bulkhead fitting tube on the water bath is connected to the pump inlet.
The water circulates but does not heat.	The temperature setting is incorrect.	Ensure that the correct set temperature has been input.
	The heater relay is defective.	Contact the Agilent Service Department for assistance. See "Obtaining Warranty and Other Services" on page 47..
	The heater elements are degraded or faulty.	For 115 V, the resistance is between 10 and 15 ohms. For 230 V, the resistance is between 48 and 52 ohms. Contact the Agilent Service Department for assistance. See "Obtaining Warranty and Other Services" on page 47.
	The main PCB is faulty.	Contact the Agilent Service Department for assistance. See "Obtaining Warranty and Other Services" on page 47.
The temperature is not stable or is incorrect.	The thermistor is faulty.	Exchange the water bath temperature probe with a known good probe if one is available. If one is not available, see " Water Bath Temperature Probe Accuracy Test " on page 38.
	Instrument requires calibration.	Contact the Agilent Service Department for assistance. See "Obtaining Warranty and Other Services" on page 47.

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

This manual contains technical reference information about the Agilent 100 Automated Disintegration Apparatus (G7962A)

- Chapter 1 Safety
- Chapter 2 Introduction
- Chapter 3 Setting Up the 100
- Chapter 4 Operating the 100
- Chapter 5 Troubleshooting and Maintenance

www.agilent.com

© Agilent Technologies, Inc. 2020, 2022
Edition: 03/2022

Document No: SD-29001863 Rev. E

