



# Thermo Scientific Digital Drybath

## Operating Manual

0000848 • Revision A • October 2021

IMPORTANT! Before using this product, read this entire operation manual carefully. Users should follow all of the operational guidelines contained in this manual and take all necessary safety precautions while using this product. Failure to follow these guidelines could result in potentially irreparable bodily harm and/or property damage.

Caution all internal adjustments and maintenance must be performed by qualified service personnel.

Material in this manual is for information purposes only. Thermo Fisher Scientific is committed to a continuing program of product development and improvement, and reserves the right to change information, such as specifications, appearance, and dimensions, described in this document without notice. Thermo Fisher Scientific makes no representations or warranties with respect to this manual. In no event shall Thermo Fisher Scientific be held liable for any damages, direct or incidental, arising out of or related to the use of this manual.

No part of this manual may be reproduced or transmitted in any form or by any means, including photocopying, recording, or using information storage and retrieval systems, for any purpose other than the purchaser's own use, without the express written permission of the manufacturer.

Any other product names and services identified in this manual are trademarks or registered trademarks of their respective owners. No such use, or the use of any trade name, is intended to convey endorsement or other affiliation with Thermo Fisher Scientific.

©2021 Thermo Fisher Scientific. All rights reserved.

All rights reserved. Without prior written consent from Thermo Fisher Scientific, any part of the manual should not be duplicated, reproduced or translated into any other languages.

Caution: All rights reserved. Thermo Fisher Scientific reserves the right to modify this manual at any time without notice.

**Before the initial use of this instrument, please carefully read this manual.**

# Contents

Models.....	1
Safety Information.....	2
Introduction .....	4
Features.....	5
Normal Working Conditions.....	5
Transportation and Storage Conditions .....	5
Basic Parameters .....	5
Regulatory Compliance .....	6
Preparation Work.....	8
Structure Diagram .....	8
Display Panel Diagram .....	8
Instructions for Buttons .....	8
Operation Guidance.....	9
Examination Before Startup.....	9
Startup .....	9
Setting Temperature and Heating Time .....	9
Start and Stop Instrument .....	10
Temperature Measurement and Calibration.....	10
Maintenance .....	11
Failure Analysis and Troubleshooting .....	12
Modular Block Accessories.....	13
Warranty .....	15
Frequently Asked Questions .....	16

# Models

This manual covers the models shown below.

**Table 1. Models**

<b>NA Model</b>	<b>EU Model</b>	<b>Voltage</b>	<b>Description</b>
88870001	N/A	100-120 Vac	TS Drybath Stdrd 1 blk 100-120 V
88870002	N/A	100-120 Vac	TS Drybath Stdrd 2 blk 100-120 V
88870003	N/A	100-120 Vac	TS Drybath Stdrd 4 blk 100-120 V
N/A	88870004	200-240 Vac	TS Drybath Stdrd 1 blk 200-240 V
N/A	88870005	200-240 Vac	TS Drybath Stdrd 2 blk 200-240 V
N/A	88870006	200-240 Vac	TS Drybath Stdrd 4 blk 200-240 V

# Safety Information

## Alert Signals



**CAUTION:** Important information is contained in any item and should be carefully read. Failure to observe the instruction would result in damage or abnormal function of the instrument.



**WARNING:** The warning message requires extremely careful operation of a certain step. Failure to observe the instruction may result in serious personal injury.

During operation, maintenance and repair of this instrument, the following basic safety notes should be observed. In case of failure to follow these instructions, the warnings or notes indicated herein, the basic protection provided by the instrument, its safety criteria of design and manufacture, and its predicted use range would be impaired. If the equipment is used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.



**CAUTION:** This instrument is for indoor use.

- **The ground connection**

To avoid the electric shock, the input power line must be reliably grounded. The instrument is equipped with the three-pin plug that has the third pin (the pin connects the ground), therefore, the plug should be used with the grounded power socket only. This is a kind of safety device. If the plug cannot be plugged into the power socket, please ask the electrician to install a correct power socket, so as to make the grounded plug to work for safety.

Thermo Fisher Scientific is not responsible for any injury as a result of the user's failure to observe the following requirements.

- **Observe voltage/power requirements**

Before the instrument is connected to the AC power source, the voltage of the power source should be the same with the required voltage of the instrument (a deviation of  $\pm 10\%$  is allowed). The rated load of the power socket should not be lower than the requirements of the instrument.

- **The power cord**

The instrument should use the power line cord attached to it. If the power line is destroyed, it must be replaced but not be repaired. The replacement should be carried out with the power line of the same type and same specification (reach out to technical service team for further information). No items should be put on the power cord when the instrument is in operation.

- **Connecting and disconnecting to/from power source**

The user should hold the plug to remove from power source. When connecting the plug, user should make sure it has been fully plugged in; when removing the plug, do not pull the power line forcefully. Do not position the equipment so that it is difficult to operate the disconnecting device (plug).

- **Placement of the instrument**

This instrument should be fixed in a low RH ( $\leq 80\%$ ) and low dust place away from water source (e.g. sink or water pipeline) and the room should be well ventilated, and free of corrosive gas or interference of strong magnetic field. The instrument should not be placed in a wet or dusty location.

The openings on this instrument are for ventilation circulation and in order to avoid over-heat of the instrument body they should not be blocked or covered. When a single set of instrument is used, the interval between ventilation opening before and after the instrument and its nearest object should not be less than 25 cm. Also, don't use the instrument on loose or soft surface, or the air inlet of instrument bottom might be blocked.

Excessive temperature will impair performance or result in failure of the instrument. This instrument should not be used in locations subjected to direct sun light. The instrument should be kept away from hot gas, oven and all other heat sources. If the instrument is to be stored for a long time, the power plug should be withdrawn and the instrument covered with soft cloth or plastic film to avoid entrance of dust. The product is powered by connecting the mains plug to a standard socket-outlet. Always place the product in such a way that it is easy for the operator to disconnect the product from the mains supply.

- **Explanation of symbols**

Attention, read user manual before use.



**WARNING:** There is a sign of “CAUTION: HOT SURFACE! ATTENTION: SURFACE HOT” on the instrument. The metal part (module) near the sign should not be touched with any part of the body when the instrument is operating in a high temperature state or just finished operation to avoid burns!



Alternating Current



Protective Conductor Terminal



**CAUTION:** In any of the following cases, immediately turn off the power supply, withdraw the power plug from the power socket, contact the supplier for service:

- Liquid drops into the inside of the instrument.
- The instrument is rained on or water is spilled on it.
- The instrument works abnormally, especially if generating an abnormal sound or odor.
- The instrument is dropped or its casing is damaged.
- The function of instrument obviously changes.

# Introduction

The product is a dry bath with advanced microcomputer control, which can be widely applied to sample reservation, enzyme reservation and reaction, DNA amplification, electrophoresis degeneration and serum coagulation, etc.

The characteristics of the product are as follows:

- The digital display for control of the temperature.
- Heat blocks help avoid possible contamination while containing the sample inside a tube.
- The heat blocks are easy to replace, clean and disinfect and are suitable for various tubes sizes.
- Built-in over-temperature protection device warrants sample and user safety.
- Temperature deviation adjustment.

# Features

This chapter introduces the usage, transportation, storage conditions of the instrument, as well as its basic parameters, performance and functions.



**CAUTION:** Before using the instrument, please make sure the working condition meets the above requirements. Especially that the power line is reliably grounded.

## Normal Working Conditions

Ambient temperature	10°C ~ 30°C
Relative humidity	≤ 80%
Power supply	100-120 V ~ 50/60 Hz or 200-240 V ~ 50/60 Hz
Altitude	Up to 2000 m
Pollution Degree	2
Overvoltage category	II Indoor use
Mains supply voltage fluctuations	±10%

## Transportation and Storage Conditions

Ambient temperature:	-20°C~+55°C
Relative humidity:	≤ 80%

## Basic Parameters

**Table 2. 100-120 V**

NA Catalog Number	88870001	88870002	88870003
Model Name	Digital 1 block 100-120 V	Digital 2 block 100-120 V	Digital 4 block 100-120 V
Temperature Range	Ambient Temperature +5°C ~ 130°C (Ambient Temperature 25°C)		
Temperature Uniformity	±1.0°C@37°C		
Temperature Accuracy	±0.5°C@37°C		
Temperature Fluctuation	±0.5°C		
Temperature Rise Time	≤ 20 min (rise from 30°C to 130°C)		≤ 25 min (rise from 30°C to 130°C)
Size (L × W × H)	288 × 200 × 100 mm 11.3 × 7.87 × 3.94 in	318 × 200 × 100 mm 12.5 × 7.87 × 3.94 in	450 × 200 × 100 mm 17.7 × 7.87 × 3.94 in
Weight	2.9 kg (6.39 lbs)	3.3 kg (7.28 lbs)	4.7 kg (10.36 lbs)
Fuse Protector (ø5×20)	F 250 V 2.5 A	F 250 V 5 A	F 250 V 8 A

**Table 3. 200-240 V**

<b>EU Catalog Number</b>	<b>88870004</b>	<b>88870005</b>	<b>88870006</b>
<b>Model Name</b>	<b>Digital 1 block 200-240 V</b>	<b>Digital 2 block 200-240 V</b>	<b>Digital 4 block 200-240 V</b>
Temperature Range	Ambient Temperature +5°C ~ 130°C (Ambient Temperature 25°C)		
Temperature Uniformity	±1.0°C		
Temperature Accuracy	±0.5°C@37°C		
Temperature Fluctuation	±0.5°C		
Temperature Rise Time	≤ 20 min (rise from 30°C to 130°C)		≤ 25 min (rise from 30°C to 130°C)
Size (L × W × H)	288 × 200 × 100 mm 11.3 × 7.87 × 3.94 in	318 × 200 × 100 mm 12.5 × 7.87 × 3.94 in	450 × 200 × 100 mm 17.7 × 7.87 × 3.94 in
Weight	2.9 kg (6.39 lbs)	3.3 kg (7.28 lbs)	4.7 kg (10.36 lbs)
Fuse Protector (ø5×20)	F 250 V 2.5 A	F 250 V 2.5 A	F 250 V 3.15 A



**CAUTION:** The instrument is only used for heating operation: from low temperature to high temperature.

## Regulatory Compliance

### European Union



The European voltage models of this product meet all the applicable requirements of the European Directives and therefore display the CE Marking. These Directives are captured in the EU Declaration of Conformity which may be obtained from the manufacturer.

### Product Safety

This product family has been tested to applicable product safety standards by a Nationally Recognized Test Laboratory (NRTL) and may bear the NRTL's mark of safety compliance to those applicable standards.

CEEMC EN 61326-1

CE Safety EN 61010-1 CE  
 Safety EN 61010-2-010  
 UL 61010-1/CSAC22.2 NO.61010-1  
 ROHS2.0 2011/65/EU

## Korean Registration



사용자안내문  
 이기기는 업무용 환경에서 사용할 목적으로 적합성 평가를 받은 기기로서 가정용 환경에서 사용하는 경우 전파 간섭의 우려가 있습니다



**WARNING:** EMC Registration is done on this equipment for business use only. It may cause interference when the product would be used in home. This warning statement applies a product for business use.

# Electromagnetic Compatibility

- Have a certified technician remove the Refractory Ceramic insulation from the unit then dispose properly.

## FCC Statement (USA)



This device complies with Part 15 Subpart B of the FCC Rules. Operation is subject to the following two conditions:

- (1) this device may not cause harmful interference, and
- (2) this device must accept any interference received, including interference that may cause undesired operation.

Any changes or modifications not expressly approved by the manufacturer could void the user's authority to operate the equipment.

**Note:** This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

## Canadian ISED IC Notice

- This ISM digital apparatus complies with Canadian ICES-001, Class A
- Cet appareil ISM est conforme à la norme NMB-001 du Canada, Classe A

## End of Life Care

Some considerations and suggestions are listed below for proper disposal of this product. While addressing these actions for safe recycling and disposal, please follow all guidelines, Safety Data Sheets (SDS), or regulations applicable to your country and region.

- This product has materials and components that may be recycled or reused according to local guidelines and regulations.
- Remove any batteries present before disposal. Batteries, battery packs, and accumulators should not be disposed of as unsorted household waste. Please use the public collection system to return, recycle, or treat them in compliance with the local regulations.
- Remove all samples and items before defrosting a unit to room ambient temperatures.
- Clean up any chemical or biological safety hazards using appropriate methods.

# Preparation Work

This chapter introduces the structure of the instrument, user interface and functions of all buttons and preparation before startup. Read the content in the chapter carefully before using the instrument for the first time.

## Structure Diagram



Figure 1. Structure Diagram

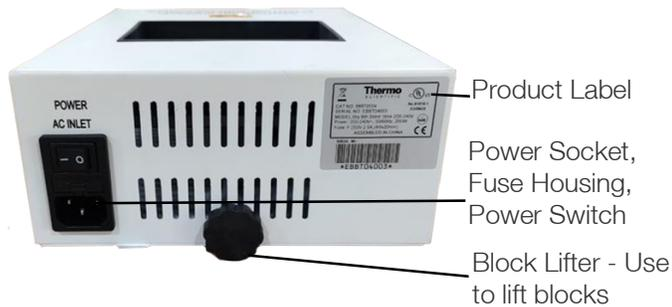


Figure 2. Back Unit

## Display Panel Diagram



Figure 3. Display Panel

## Instructions for Buttons

	<p>Setting or adjusting button: Press the button to display present setting temperature and press the button again to display present time. After the unit has reached set point and temperature is stable, calibration can be set by pressing the button.</p>
	<p>Under the status of setting, press “▲” or “▼” to modify the flashing figure. The instrument will enter temperature calibration mode if user press the two buttons simultaneously under the status of stopping.</p>
	<p>When setting or adjusting, pressing the button can move the flashing figure. During the operation, the display can be switched between the temperature and the time.</p>
	<p>For starting up or stopping program.</p>

# Operation Guidance

This chapter introduces the displaying window and the operation of buttons of the dry bath with constant temperature.

**Note:** The illumination pattern of the 'Heat Light' indicates the Operation status.

- Heat Light OFF (Unit Stopped) = "Set/Cal" button used for Setpoint adjustment.
- Heat Light blinks/flashes as Block is heating or cooling toward setpoint.
- Heat Light ON (Unit Started and stable at setpoint) = "Set/Cal" used for Calibration.

## Examination Before Startup

Please confirm the following content before plugging the plug into the power socket.

1. The power source is in accordance with the voltage requirement of the instrument (please refer to the chapter II for the requirement of the voltage).
2. Make sure the plug has been fully plugged into the power socket.
3. The power line is reliably grounded.



**CAUTION:** If the display of the instrument is abnormal after startup, please turn off the power source immediately and contact with the supplier.

## Startup

Press the power switch and the instrument will power on. All indicator lights and display serial ports will be turned on. About 3 ~ 5 seconds later, temperature display window will show the real-time temperature. All indicator lights will be turned off.

## Setting Temperature and Heating Time

Set Temperature and Hold-Time Setpoint:

1. Connect power cord to appropriate electrical supply.
2. Make Power switch, located near power cord on rear, to the 'ON' position.
3. Confirm the Heat-indicator is 'off'. If 'ON' or 'blinking', press "Stop".
4. Set-up Temperature:
  - a. Press the "SET/CAL" button. The instrument is in the Temperature Setpoint mode with the display window showing the previous temperature set value, and the left digit flashing.
  - b. Press "▲" to increase or "▼" to decrease the flashing digit of the Setpoint.
  - c. Press "SHIFT" to move the flashing right, to the next digit.
  - d. Repeat Steps 4b and 4c to complete Temperature setpoint selection.
5. Setting Hold-Time Setpoint:
  - a. After the temperature is selected, press "SET/CAL" button. The instrument is in the Hold-time Setpoint mode with the display window showing the previous Hold-time set value, and the left digit flashing.
  - b. Press "▲" to increase or "▼" to decrease the flashing digit of the Hold-Time. Choose a time from 00:01 to 99:59, hours and minutes.
  - c. Press "SHIFT" to move the flashing right, to the next digit.
  - d. Repeat Steps 5b and 5c to complete Hold-Time setpoint selection. For Continuous operation, Choose 99:59 then increase to display "- :- -".
  - e. After the Time setting is complete, Wait for 5 seconds and the instrument will automatically exit the Setpoint mode. The display will now show the current measured temperature.
6. Press "Start" button to begin heating the Block. The green "Heat Light" will blink as Block is heating toward set point.
7. Allow 30-minutes for temperature to stabilize. The heat indicator stops blinking and remains constantly "ON" to indicate the temperature is stable, and instrument is ready to use.

**Instruction:** There is limitation for temperature and time setting. The setting temperature range is: 15 ~ 135°C; the time range is: 0 ~ 99 h 59 min and continuous mode "- -:- -".

## Start and Stop Instrument

**START:** Press "START/STOP" to start operation, the indicator light "HEAT" will be flashing. When the instrument achieves the setting temperature, the "HEAT" light will stop flashing. This indicates that the unit will enter hold temperature status. The digital display window will display the real-time temperature and the timing function by pressing "SHIFT" button. When the holding time equals to the setting time, the instrument will alarm 5 beeps.

**STOP:** Under the working status, press "START/STOP" to stop working. At that time, corresponding working indicator light "HEAT" will be turned off. The digital display window displays the real-time temperature.

## Temperature Measurement and Calibration

1. The instrument has been operational for minimum of 30-minutes at desired setpoint.
2. An independent temperature measurement device is in the designed 5 mm hole in Block, along with paraffin oil, or similar heat-transfer oil.
3. Confirm the Heat-indicator is 'ON' steady, not 'blinking'.
4. Compare the Display and independent measurements. And check if the temperature gap is  $> 0.2^{\circ}\text{C}$  or not.
5. If less than  $0.2^{\circ}\text{C}$ , no calibration should be made.
6. If more than  $0.2^{\circ}\text{C}$ , calibrate the Temperature by these steps:
  - a. Press and hold the "SET/CAL" button. The instrument is in the Temperature Calibration mode with the display window showing the measured temperature value, and the left digit flashing.
  - b. Press "▲" to increase, or "▼" to decrease, the flashing digit to match the same digit of the independent measurement.
  - c. Press "SHIFT" to move the flashing right, to the next digit.
  - d. Repeat Steps 6b and 6c to complete Temperature calibration.
  - e. After the temperature calibration is completed, press "SET/CAL" button to exit the calibration mode and continue the Start mode.

7. The Heat-indicator may start 'blinking' as the temperature is stabilized to the calibrated temperature.
8. When the Heat-indicator is 'ON' steady, the temperature is stable.
9. After the instrument stable operated for additional 10 minutes, compare the display temperature and measurement again.  
If the temperature gap is  $> 0.2^{\circ}\text{C}$ , repeat all steps of 6.



### CAUTION:

- To ensure the accuracy of temperature, the instrument shall be calibrated after temperature stabilizes for 30 minutes.
- Calibrate the instrument with qualified Grade II standard filled thermometer.
- Calibration point: Fill the middle hole of the heat block with paraffin oil and soak the thermometer bulb.

# Maintenance

The holes of the blocks should be regularly cleaned with the damp cloth to ensure the test tube be well contacted with the wall of the holes so as to have good heat conduction. If the surface of the instrument is polluted, it can be cleaned with a slightly damp soft cloth. Generally unit falls under protective system of IP20 pursuant to IEC 60529.



**WARNING:** When cleaning the instrument, the power supply should be shut off and unplugged. The instrument surface should be cleaned with a non-corrosive cleaning agent.

# Failure Analysis and Troubleshooting

This chapter introduces possible failures of this instrument, explanation and troubleshooting.



**CAUTION:** The user is not allowed to open the housing of the instrument for inspection during the warranty period. If any failures that require opening the housing for inspection occur, the supplier or manufacturer should be contacted timely.

**Table 4. Failure Analysis and Troubleshooting**

No.	Error	Cause	Solution
1	Display window is not turned on after switching on power supply	Power supply fails to be connected	Check power supply and connect it.
		Fuse burned out	Replace fuse
		Damage of the switch	Replace the switch.
		Others	Contact Thermo Fisher Scientific.
2	Temperature display window shows "001" and the buzzer alarm in a di...di...sound	Short circuit of sensor	Check if the connecting line of sensor is damaged, and replace it.
3	Temperature display window shows "002" and the buzzer alarm in a di...di...sound	Open circuit of sensor	Check if the connecting line of sensor is damaged, and replace it.
4	Temperature display does not conform to actual temperature	Sensor is damaged or in poor contact	Contact Thermo Fisher Scientific.

# Modular Block Accessories

Table 5. Modular Block Accessories

	Catalog No.	Heating Blocks	No. of Holes	Block Dimensions W x D x H (mm/in)
	88870101	For 6 mm test tubes	46	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870102	For 10 mm test tubes	28	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870103	For 1.5 ml test tubes	28	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870104	For 2.0 ml test tubes	28	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870105	For 12 mm & 13 mm test tubes	24	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870106	For 15 mm & 16 mm test tubes	15	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870107	For 17 mm & 18 mm test tubes	12	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870108	For 20 mm test tubes	8	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870109	For 25 mm test tubes	6	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870110	For 0.5 ml test tubes	40	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870111	For 96 well Elisa plate	1	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870112	For mixed size test tubes 32 holes for 6 mm test tubes 21 holes for 10 mm test tubes	32 + 21	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870113	For mixed size test tubes 18 holes for 1.5 ml test tubes 10 holes for 2.0 ml test tubes	18 + 10	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870114	For mixed size test tubes 3 holes for 25 mm test tubes 12 holes for 13 mm test tubes 6 holes for 6 mm test tubes	3+12+6	124 x 76 x 39 / 4.9 x 3.0 x 1.5

**Table 5. Modular Block Accessories**

	<b>Catalog No.</b>	<b>Heating Blocks</b>	<b>No. of Holes</b>	<b>Block Dimensions W x D x H (mm/in)</b>
	88870115	For mixed size test tubes 30 holes for 0.5 ml test tubes 20 holes for 0.2 ml test tubes	30+20	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870116	For 15 ml flat bottom test tubes	15	124 x 76 x 113 / 4.9 x 3.0 x 4.55
	88870117	For 50 ml flat bottom test tubes	4	124 x 76 x 113 / 4.9 x 3.0 x 4.55
	88870118	For 15 ml conical bottom test tubes	15	124 x 76 x 113 / 4.9 x 3.0 x 4.55
	88870119	For 50 ml conical bottom test tubes	4	124 x 76 x 113 / 4.9 x 3.0 x 4.55
	88870120	For 96 well Non-skirted PCR plate	1	124 x 76 x 39 / 4.9 x 3.0 x 1.5
	88870121	For 96 well half/full skirted PCR plate	1	124 x 76 x 55 / 4.9 x 3.0 x 2.17
	88870122	Temperature probe, PT1000, stainless steel	NA	NA
* Dry bath blocks can be autoclaved.				

# Warranty

When used in laboratory conditions and according to these operation instructions and maintenance, this product is warranted for 24 months against defective materials or workmanship. The 24 month warranty period begins from the delivery date of this product.

For product quality or performance issues, contact Thermo Fisher Scientific Customer Service.

# Frequently Asked Questions

## Scenario 1

Does unit turn off once the program time is complete:

Example- If unit is set for one program to run at 45°C for 2 hrs.

**Q**-Would the unit turn off upon completion of this 2 hrs?

**A**-The instrument doesn't turn off by itself, it just stops running.

**Q**-If not- would the unit hold the set temperature beyond this 2 hrs?

**A**-After this 2 hrs, the unit doesn't hold the set temperature. It will cool to ambient temp.

**Q**-Would unit return to ambient temp?

**A**-The unit will return to ambient temp.

## Scenario 2

Suppose the unit is running a set program and there is power outage. Once the power resumes-

**Q**-Would unit restart automatically once power is restored

**A**-The instrument will automatically turn on.

**Q**-Would unit resume the program from where it was left?

**A**-It can't.

**Q**-Would customer need to start the program manually?

**A**-Yes.

**Q**-Would the unit keep the saved program in memory?

**A**-The unit has no memory function.

**Q**-Would customer need to create new program?

**A**-Need to reset the program.

# WEEE Compliance

**WEEE Compliance.** This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Thermo Fisher Scientific products which may assist the detection of substances subject to the RoHS Directive are available at [thermofisher.com/WEEERoHS](http://thermofisher.com/WEEERoHS).

## Great Britain



**WEEE Konformität.** Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliedstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, dss3A die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter [thermofisher.com/WEEERoHS](http://thermofisher.com/WEEERoHS).

## Deutschland



**Conformità WEEE.** Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2012/19/EU in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo Fisher Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito [thermofisher.com/WEEERoHS](http://thermofisher.com/WEEERoHS).

## Italia



**Conformité WEEE.** Ce produit doit être conforme à la directive européenne (2012/19/EU) des Déchets d'Équipements Électriques et Électroniques (DEEE). Il est marqué par le symbole suivant. Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur [thermofisher.com/WEEERoHS](http://thermofisher.com/WEEERoHS).

## France





### **North America's Technical Support Team**

Phone: 1-866-984-3766, Option#2

Service Email: [Servicesupport.led.asheville@thermofisher.com](mailto:Servicesupport.led.asheville@thermofisher.com)

Presales Email: [LED.Presales@thermofisher.com](mailto:LED.Presales@thermofisher.com)

Webpage: [www.unitylabservices.com](http://www.unitylabservices.com)

### **Europe Middle East Africa Technical Support Team**

German national toll free: 0800 1 112 110, Fax: 0800 1 112 114

German international: +49 6103 408 0, Fax: +49 6103 408 12 13

Email: [unity.de@thermofisher.com](mailto:unity.de@thermofisher.com) and [let.support@thermofisher.com](mailto:let.support@thermofisher.com)

### **Singapore Technical Support Team**

Service Hotmail is [Service.LPG.ROA1@thermofisher.com](mailto:Service.LPG.ROA1@thermofisher.com)

Thermo Fisher Scientific Inc.  
22 Alpha Rd,  
Chelmsford, MA  
01824, USA

Find out more at [thermofisher.com/drybaths](http://thermofisher.com/drybaths)

Australia: (613) 9757-4300 In Australia: (1300) 735-295  
China: (86) 21-6865-4588 Germany: (49) 6184-90-6321  
India: (91) 22-4157-8800 Japan: (81) 045-453-9175  
North America: 1-978-232-6000 Toll Free: 1-800-225-1480  
Singapore: (65) 6778-6876



*Thermo Fisher Scientific  
Water Analysis Instruments  
Chelmsford, MA USA  
Quality Management System  
Registered to ISO 9001*

© 2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. 1021

**ThermoFisher**  
S C I E N T I F I C