



Thermo Scientific Touch Screen Drybath

Operating Manual

0000849 • 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

Preface	1
Models.....	2
Safety Information.....	3
Introduction	5
Features.....	6
Normal Working Conditions.....	6
Transportation and Storage Conditions	6
Basic Parameters	6
Regulatory Compliance	7
Preparation Work.....	9
Structure Interface.....	9
Operation Guidance.....	10
Examination Before Start-Up	10
StartUp	10
Operating Instructions	10
Maintenance	14
Failure Analysis and Troubleshooting	15
Modular Block Accessories.....	16
Warranty	18
Warranty Scope	18
FAQ Section	19

Preface

Our **Sales Support** staff can provide information on pricing and give you quotations. We can take your order and provide delivery information on major equipment items or make arrangements to have your local sales representative contact you. Our products are listed on the Internet and we can be contacted through our Internet home page.

Our **Service Support** staff can supply technical information about proper setup, operation or troubleshooting of your equipment. We can fill your needs for spare or replacement parts or provide you with on-site service. We can also provide you with a quotation on our Extended Warranty for your Thermo Scientific products.

Whatever Thermo Scientific products you need or use, we will be happy to discuss your applications. If you are experiencing technical problems, working together, we will help you locate the problem and, chances are, correct it yourself over the telephone without a service call.

When more extensive service is necessary, we will assist you with direct factory trained technicians or a qualified service organization for on-the-spot repair. If your service need is covered by the warranty, we will arrange for the unit to be repaired at our expense and to your satisfaction.

Regardless of your needs, our professional telephone technicians are available to assist you Monday through Friday from 8:00 a.m. to 6:00 p.m. Eastern Time. Please contact us by telephone or fax.

International customers, please contact your local Thermo Scientific distributor.

Models

This manual covers the models shown below.

Table 1. Models

NA Model	EU Model	Voltage	Description
88870007	N/A	100-120 Vac	Drybath Stdrd 1 blk 100-120 V
88870008	N/A	100-120 Vac	Drybath Stdrd 2 blk 100-120 V
88870009	N/A	100-120 Vac	Drybath Stdrd 4 blk 100-120 V
N/A	88870010	200-240 Vac	Drybath Stdrd 1 blk 200-240 V
N/A	88870011	200-240 Vac	Drybath Stdrd 2 blk 200-240 V
N/A	88870012	200-240 Vac	Drybath Stdrd 4 blk 200-240 V

Safety Information

Alert Signals



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



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

Note: Notes alert you to pertinent facts and conditions.



WARNING: If the Thermo Scientific Touch Screen Drybath is used in a manner not specified by the manufacturer, the protection provided by the unit may be impaired.

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 here in, 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.

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

For your own safety, user should observe the following safety warning signs.

The warning signs indicate potential sources of danger. They also inform on how correct action can avert danger.



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.

- **Observe voltage/power requirements**

Before the instrument is connected with the AC power source, the voltage of the power source should be the same with the required voltage of the instrument (a deviation of 10% is allowed). e 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 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 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 instrument. This instrument should not be used in location 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!



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.



Alternating Current

Introduction



CAUTION: After unpacking, immediately check the goods against the packing list. In case of damaged or missed goods, please immediately contact the distributor in which the product was purchased from. After unpacking, the packing box and packing materials should be well kept for future repair. For the damage of instrument occurring during transporting to the repair department due to improper package, Thermo Fisher Scientific will not be held responsible for damages.

The product is dry bath with advanced microprocessor 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 and control of the temperature.
- Heat blocks help avoid possible contamination while containing the sample inside a tube.
- The heat blocks are easily replaced, cleaned or disinfected and are suitable for various tubes.
- 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, make sure the working conditions meet the above requirements.

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

Parameters / Type	Touch Screen 1 block 100-120 V	Touch Screen 2 block 100-120 V	Touch Screen 4 block 100-120 V
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)	373 × 200 × 100 mm 4.7 × 7.87 × 3.94 in	403 × 200 × 100 mm 15.9 × 7.87 × 3.94 in	535 × 200 × 100 mm 21.1 × 7.87 × 3.94 in
Weight	3.4 kg (7.5 lbs)	3.7 kg (8.16 lbs)	5.1 kg (11.24 lbs)
Fuse Protector (ø5×20)	F 250 V 2.5 A	F 250 V 5 A	F 250 V 8 A

Table 3. 220 V

Parameters / Type	Touch Screen 1 block 200-240 V	Touch Screen 2 block 200-240 V	Touch Screen 4 block 200-240 V
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)	373 × 200 × 100 mm 14.7 × 7.87 × 3.94 in	403 × 200 × 100 mm 15.9 × 7.87 × 3.94 in	535 × 200 × 100 mm 21.1 × 7.87 × 3.94 in
Weight	3.4 kg (7.5 lbs)	3.7 kg (8.16 lbs)	5.1 kg (11.24 lbs)
Fuse Protector (ø5×20)	F 250 V 2.5 A	F 250 V 2.5 A	F 250 V 3.15 A

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.

Korean Registration



사용자안내문
이기는은무용환경에서 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.

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

Electromagnetic Compatibility

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

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 startup when using the instrument for the first time.

Structure Interface

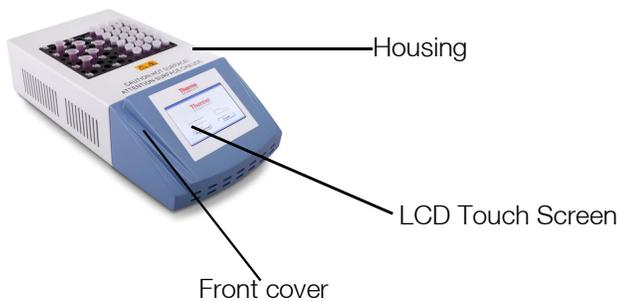


Figure 1. Unit Front

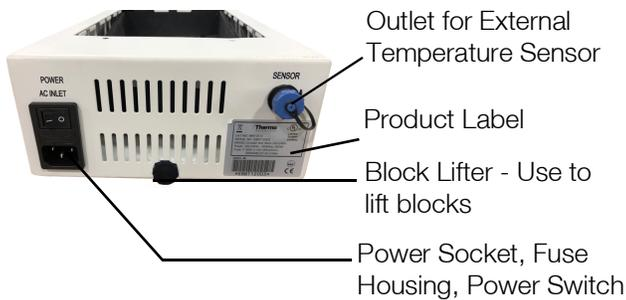


Figure 2. Unit Back

Note:

- When the external temperature sensor is connected, the internal temperature sensor will disconnect automatically, and the temperature control and display will relay on the external temperature sensor.
- DO NOT plug in or remove the external temperature sensor without the power being off.

Operation Guidance

The chapter introduces the display window and the operation of buttons of the dry bath with constant temperature.

Examination Before Start-Up

Confirm the following before inserting the plug into the power socket.

1. Ensure the power source is in accordance with the voltage requirement of the instrument (refer to “Features” section for voltage requirements).
2. Make sure the plug has been completely inserted into the power socket.
3. Ensure the power line has been reliably grounded.



CAUTION: If the display of the instrument is abnormal after startup, turn off the power source immediately and contact Thermo Fisher Scientific.

StartUp

Press the power switch and the instrument will be powered on. The LCD screen of the operation panel will show the main interface.

Operating Instructions

The Main Interface

- Press “Single point set” to enter into the single setting interface.

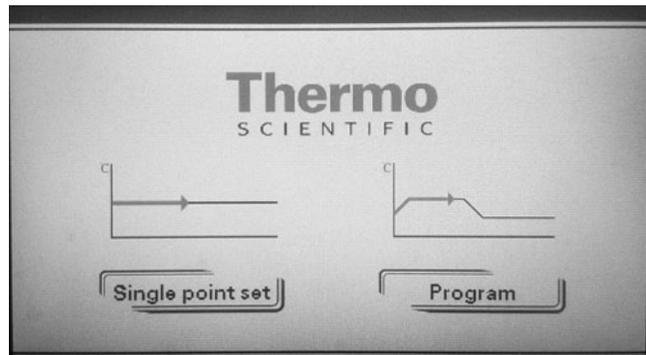


Figure 3. Main Interface

Single Setting Interface

- Press “BACK” to go back to the main interface.
- Press “RUN” to enter into the single running interface.
- Press “Temp” to set the temperature.
- Press “Time” to set the time.



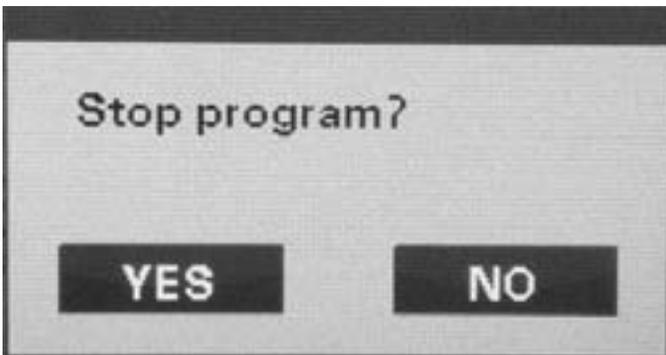
Figure 4. Single Setting Interface

Instruction: There is limitation for temperature and time setting. The setting temperature range is: 15~135°C; the time range is: 0~99: 59 and “, ”.

Single Running Interface



- Press “STOP”, and the interface of asking you whether to stop the program will appear.



- Press “YES” to stop the program and to go back to the main interface.
- Press “NO” to go back to the single running interface.

Calibration

- Press “CAL” to calibrate the temperature adjustment interface. (If the external sensor is being used, then there would be no “CAL.”, and the temperature is unable to be adjusted.)



Figure 5. Calibrating the temperature

- Input the correct temperature into the box that pops out.
- To ensure the accuracy of temperature, the instrument should be calibrated after temperature stabilizes for 30 minutes.
- Calibrate the instrument with the qualified Grade II standard filled thermometer.
- Calibration point: The middle hole of the heat block. Pour paraffin oil into the well, and soak the thermometer bulb.

Instruction: During the operation of the instrument, the temperature and the time should not be changed. If it they should, stop the program and enter into the setting interface to change them.

Listed Files Interface

- Press “BACK” to go back to the main interface.
- Enter into the setting interface of the specific file in accordance with its name. If the file has a password, then enter into the password inputting interface.

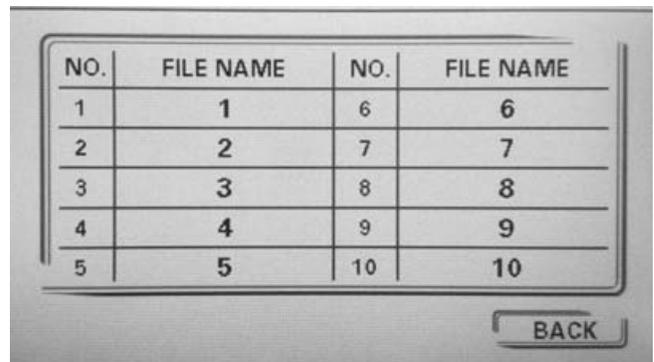


Figure 6. Listed Files Interface

Password Inputting Interface

- Input the password into the box.



CAUTION: If the password has been forgotten, user can use 000000 to ROOT and then the file will be reset to the initial state (the name of the file would not be changed).



Figure 7. Password Interface

Program Setting Interface

- Press “BACK” to go back to the listed files interface.
- Press “SAVE” to enter into the files saving interface
- Press “RUN” to enter into the program running interface. If the alteration of the program has not been saved yet, the warning box will pop up.



Figure 8. Program Setting Interface



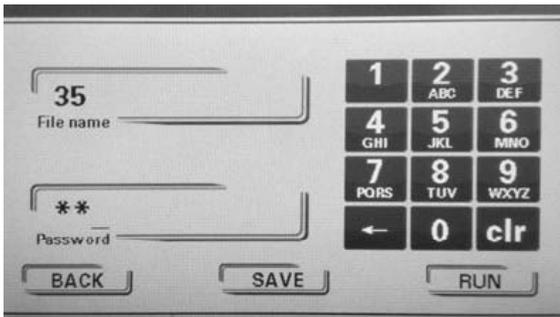
- Press “YES”, and the program will run in the changed form, but the changes won’t be saved. Press “NO” to go back to the program setting interface.
- The user can set 1-5 sections of the program. Press  to delete the corresponding section (when there is only one section, this step would be of no use).
- Press  to add sections. Choose the corresponding box to set the corresponding temperature and time.



CAUTION: This instrument doesn't have cooling function. If the user wants to set multiple sections, set the temperature from low to high. Otherwise, the temperature will be very slowly decreasing while the blocks cool down.

Files saving interface

- Press “BACK” to go back to the program setting interface.
- Press “SAVE” to save the file.
- Press “RUN” to enter into the program running interface. If the changed program has not been saved, the warning box will pop up.
- Choose “File name” and “Password” to set the file name and the password. By pressing Figures from 2-9 quickly, the input method will be changed to English to set the file name.



Instruction: During the operation of the instrument, the temperature and the time should not be changed. If needed, stop the program and enter into the setting interface to change them.

Figure 9. Files Saving Interface

Program Running Interface

- Press “CAL.” to call for the temperature adjustment interface. (If the external sensor is being used, then there would be no “CAL”, and the temperature is unable to be adjusted). Refer to ‘Single Running Interface’.
- Press “SKIP” to skip to the next section. If it is the last section, the interface of asking you whether to stop the program will appear. Refer to ‘Single Running Interface’.



Figure 10. Program Running Interface

- Press “PAUSE”, the program will be paused and the temperature will be maintained to the set temperature and the pause interface will pop up.

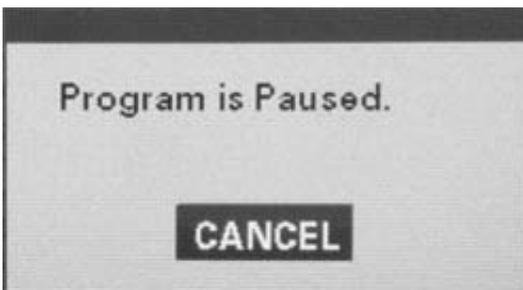


Figure 11. Program Paused

- Press “STOP”, the cancel interface will pop up. Refer to ‘Single Running Interface’.

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 details possible failures of this instrument and possible causes with troubleshooting.

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	Replace fuse
		Damage of the switch	Replace the switch
		Others	Contact Thermo Fisher Scientific
2	The LCD shows “Err code: 0x1” and the buzzer alarms in a di...di...sound	Short circuit of sensor	Check whether the connecting wire of sensor is damaged and replace it
3	The LCD shows “Err code: 0x2” and the buzzer alarms in a di...di...sound	Open circuit of sensor	Check whether the connecting wire of sensor is damaged, and replace it
4	The LCD shows “Err code: 0x3” and the buzzer alarms in a di...di...sound	Sensor is damaged or the heat blocker is damaged	Contact supplier or factory
		Use the external sensor without putting it into the module	Put the external sensor correctly into the corresponding place of the module
5	Temperature display does not conform to actual temperature	Sensor is damaged or in poor contact	Contact supplier or factory



CAUTION: The user is not allowed to open the housing of the instrument for inspection during the warranty period. If any failures result from opening the housing for inspection, 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

Table 5. Modular Block Accessories

	Catalog No.	Heating Blocks	No. of Holes	Block Dimensions W x D x H (mm/in)
	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
	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 none 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	N/A	NA
* Dry bath blocks can be autoclaved.				

Warranty

The Warranty Period starts two weeks from the date your equipment is shipped from our facility. This allows for shipping time so the warranty will go into effect at approximately the same time your equipment is delivered. The warranty protection extends to any subsequent owner during the first year warranty period.

During the first two (2) years, component parts proven to be non-conforming in materials or workmanship will be repaired or replaced at Thermo's expense, labor included. Installation and calibration are not covered by this warranty agreement. The Technical Services Department must be contacted

for warranty determination and direction prior to performance of any repairs. Expendable items, glass, filters and gaskets are excluded from this warranty.

Replacement or repair of components parts or equipment under this warranty shall not extend the warranty to either the equipment or to the component part beyond the original warranty period. The Technical Services Department must give prior approval for return of any components or equipment.

At Thermo's option, all non-conforming parts must be returned to Thermo Fisher Scientific postage paid and replacement parts are shipped FOB destination.

THIS WARRANTY IS EXCLUSIVE AND IN LIEU OF ALL OTHER WARRANTIES, WHETHER WRITTEN, ORAL OR IMPLIED. NO WARRANTIES.

OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE SHALL APPLY. Thermo shall not be liable for any indirect or consequential damages including, without limitation, damages relating to lost profits or loss of products.

Your local Thermo Sales Office is ready to help with comprehensive site preparation information before your equipment arrives. Printed instruction manuals carefully detail equipment installation, operation and preventive maintenance.

Warranty Scope

The above mentioned warranty does not apply to damages incurred by improper use and maintenance, using under unqualified conditions, unauthorized maintenance or modification.

Thermo Fisher Scientific does not provide any other warranty expressions otherwise; and does not bear liabilities for business promotion under special circumstances.

FAQ Section

:Scenario 1: Does unit turn off once the program time is complete: Example- If unit is set for one program to run at 45degC 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.

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, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter 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.

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.

France





North America's Technical Support Team

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

Service Email: Servicesupport.led.asheville@thermofisher.com

Presales Email: LED.Presales@thermofisher.com

Webpage: 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 and let.support@thermofisher.com

Singapore Technical Support Team

Service Hotmail is Service.LPG.ROA1@thermofisher.com

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

Find out more at 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