



Modular Block Dri-Baths

Microprocessor Control:

2000Q, 2000-1CEQ
2001Q, 2001-1CEQ
2002Q, 2002-1CEQ
2003Q, 2003-1CEQ
2004Q, 2004-1CEQ

Bimetallic Control:

2050Q, 2050MARQ, 2050SUPQ, 2050IDQ
2050-1CEQ, 2050-1CEMARQ, 2050-1CESUPQ, 2050-1CEIDQ
2052Q, 2052IDQ, 2052-1CEQ, 2052-1CEIDQ
2053Q, 2053-1CEQ,
2054Q, 2054IDQ, 2054-1CEQ, 2054-1CEIDQ
2056Q, 2056-1CEQ

Table of Contents

Safety Information	3
Alert Signals.....	3
Introduction.....	4
Specifications	5
Digital Units	5
Analog Units	6
Declaration of Conformity	7
Unit's Environmental Operating Conditions	7
Power Requirements	8
Unpacking and Installation.....	9
Shipping Carton	9
Location	9
Heating Blocks.....	9
Electrical Requirements.....	9
Operation	10
Power Up.....	10
Position Block for Temperature Sensor Probe, 2000 Series	10
Control Panel: 2000 Series	11
Setting the Temperature: 2000 Series	11
Temperature Calibration: 2000 Series	12
Control Panel: 2050 Series	13
Setting the Temperature: 2050 Series	13
Maintenance.....	15
Troubleshooting	16
Replacement Parts.....	17
Accessories	18
Ordering Procedures	19
Warranty.....	20

Safety Information

Alert Signals

**Warning**

Warnings alert you to a possibility of personal injury.

**Caution**

Cautions alert you to a possibility of damage to the equipment.

**Note**

Notes alert you to pertinent facts and conditions.

**Hot Surface**

Hot surfaces alert you to a possibility of personal injury if you come in contact with a surface during use or for a period of time after use.

**Warning**

If the Thermo Scientific Modular Block Dri-bath is used in a manner not specified by the manufacturer, the protection provided by the unit may be impaired.

Thank you for selecting Thermo Scientific for your equipment needs.

Your Thermo Scientific Modular Block Dri-bath has been designed with function, reliability, and safety in mind. It is your responsibility to install it in conformance with local electrical codes. It is most important that the user follow installation instructions exactly as written. Do not attempt operation without this information.

Introduction

The Thermo Scientific Modular Block Dri-Baths provide uniform dry heating for cuvettes, microcentrifuge tubes, titer plates and test tubes of various sizes. Close contact of tubes to block walls provides excellent heat transfer. These dri-baths are designed for use in food sciences, chemistry, hematology, pathology or clinical labs and are ideal for general heating and denaturing.

Large selections of interchangeable heating blocks provide the dri-baths with the flexibility of accommodating various test tube sizes and quantities. In addition, a solid block permits the heaters to be used as a temperature-controlled hotplate. The cabinets are made from steel with a powder-coated finish. Heating blocks are solid aluminum with a sealed black anodized finish for superior heat absorption.

A PID microprocessor-based controller maintains temperature from slightly above ambient to 130°C for 2000 series. A patented temperature sensing probe system ensures accurate block temperature by measuring the temperature within the block.

Temperature is maintained from slightly above ambient to 130°C by 2 control thermostats for 2050 series, one for lower (warm) temperatures and another for higher (hot) temperatures. Temperature uniformity in the heating block is maintained to within $\pm 5^{\circ}\text{C}$. The 3-position rocker switch can be set LOW, HIGH or OFF to produce the required temperature range. A separate heater status lamp indicates when power is being supplied through either thermostat to the heating element.

Specifications

Digital Units

Part Number	Temperature Range	Temp. Control @37°C	Temp. Uniformity @37°C	Unit's Overall Dimensions	Heater Block Dimensions	Block Capacity	Shipping Weights
2000Q	Slightly above ambient to 130° C	± 0.5° C	±0.4°C	8-9/16" x 7-1/2" x 3-3/8" (22 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	1 block	(2 kg)
2000-1CEQ	Slightly above ambient to 130° C	±0.5° C	±0.4° C	8-9/16" x 7-1/2" x 3-3/8" (22 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	1 block	5 lbs. (2 kg)
2001Q	Slightly above ambient to 130° C	±0.5° C	±0.4° C	8-9/16" x 7-1/2" x 3-3/8" (22 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	2 block	(3 kg)
2001-1CEQ	Slightly above ambient to 130° C	0.5° C	±0.4°C	8-9/16" x 7-1/2" x 3-3/8" (22 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	2 block	6 lbs. (3 kg)
2002Q	Slightly above ambient to 130° C	0.5° C	±0.4°C	8-9/16" x 13-1/2" x 3-3/8" (22 x 34 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	3 block	(4 kg)
2002-1CEQ	Slightly above ambient to 130° C	0.5° C	±0.4°C	8-9/16" x 13-1/2" x 3-3/8" (22 x 34 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	3 block	(4 kg)
2003Q	Slightly above ambient to 130° C	0.5° C	±0.4°C	(28 x 40 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	4 block	(4 kg)
2003-1CEQ	Slightly above ambient to 130° C	0.5° C	±0.4°C	(28 x 40 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	4 block	(4 kg)
2004Q	Slightly above ambient to 130° C	0.5° C	±0.4°C	(28 x 40 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	6 block	(5 kg)
2004-1CEQ	Slightly above ambient to 130° C	0.5° C	±0.4°C	(28 x 40 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	6 block	(5 kg)

SPECIFICATIONS

Analog Units

Part Number	Temperature Range	Temp. Control @37°C	Temp. Uniformity @37°C	Unit's Overall Dimensions	Heater Block Dimensions	Block Capacity	Shipping Weights
2050Q	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	6-3/16" x 7-5/8" x 3-3/8" (16 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	1 block	5 lbs. (2 kg)
2050-1CEQ	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	6-3/16" x 7-5/8" x 3-3/8" (16 x 19 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	1 block	(2 kg)
2052Q	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	8-5/16" x 8-5/8" x 3-3/8" (21 x 22 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	2 block	6 lbs. (3 kg)
2052-1CEQ	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	8-5/16" x 8-5/8" x 3-3/8" (21 x 22 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	2 block	(3 kg)
2053Q	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	11-1/4" x 9-1/4" x 3-3/8" (29 x 24 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	3 block	(4 kg)
2053-1CEQ	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	11-1/4" x 9-1/4" x 3-3/8" (29 x 24 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	3 block	(4 kg)
2054Q	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	8-5/16" x 12-3/8" x 3-3/8" (21 x 31 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	4 block	(4 kg)
2054-1CEQ	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	8-5/16" x 12-3/8" x 3-3/8" (21 x 31 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	4 block	(4 kg)
2056Q	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	11-1/4" x 12-7/8" x 3-3/8" (29 x 33 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	6 block	(5 kg)
2056-1CEQ	Slightly above ambient to 130° C; Low range, slightly above ambient to 60° C; High range, 50° C to 130° C	±3.5°C	±0.5°C	11-1/4" x 12-7/8" x 3-3/8" (29 x 33 x 9 cm)	3-3/4" x 2-15/16" x 2" (9.5 x 7.5 x 5.1 cm)	6 block	(5 kg)

Note: High end of low range can reach 80°C.

High end of high range can reach 160°C.

Declaration of Conformity

(for 220-240 volt, -33 CE models only)

We hereby declare under our sole responsibility that this product conforms with the technical requirements of the following standards:

EMC:	EN 61000-3-2	Limits for harmonic current emissions
	EN 61000-3-3	Limits for voltage fluctuations and flicker
	EN 61326-1	Electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
Safety:	EN 61010-1	Safety requirements for electrical equipment for measurement, control, and laboratory use; Part I: General Requirements
	EN 61010-2-051	Part II: Particular requirements for laboratory equipment for the heating of materials

per the provisions of the Electromagnetic Compatibility Directive 2004/108/EC, and per the provisions of the Low Voltage Directive 2006/95/EC.

The authorized representative located within the European Community is:

Thermo Fisher Scientific
Electrothermal House
Unit 12A
Purdeys Industrial Estate
Purdeys Way
Rochford, Essex, SS4 1ND
United Kingdom
Tel: +44(0)1702 303350

Copies of the Declaration of Conformity are available upon request.

Unit's Environmental Operating Conditions

POLLUTION DEGREE*:	2
INSTALLATION CATEGORY*:	II
ALTITUDE:	2000 Meters MSL (Mean Sea Level)
HUMIDITY:	80% maximum, non-condensing
ELECTRICAL SUPPLY:	120VAC or 240VAC
VOLTAGE TOLERANCE:	±10% of normal rated line
TEMPERATURE:	15°C to 40°C
PRODUCT USAGE:	This product is intended for use indoors only

SPECIFICATIONS

Power Requirements

2000Q:	120 VAC, 50/60 Hz, 100 Watts, 0.83 Amps
2000-1CEQ:	240 VAC, 50/60 Hz, 100 Watts, 0.42 Amps
2001Q:	120 VAC, 50/60 Hz, 150 Watts, 1.25 Amps
2001-1CEQ:	240 VAC, 50/60 Hz, 150 Watts, 0.63 Amps
2002Q:	120 VAC, 50/60 Hz, 200 Watts, 1.67 Amps
2002-1CEQ:	240 VAC, 50/60 Hz, 200 Watts, 0.80 Amps
2003Q:	120 VAC, 50/60 Hz, 200 Watts, 0.83 Amps
2003-1CEQ:	240 VAC, 50/60 Hz, 300 Watts, 1.25 Amps
2004Q:	120 VAC, 50/60 Hz, 400 Watts, 3.33 Amps
2004-1CEQ:	240 VAC, 50/60 Hz, 400 Watts, 1.67 Amps
2050Q:	120 VAC, 50/60 Hz, 100 Watts, 0.83 Amps
2050MARQ:	120 VAC, 50/60 Hz, 100 Watts, 0.83 Amps
2050SUPQ:	120 VAC, 50/60 Hz, 100 Watts, 0.83 Amps
2050IDQ:	120 VAC, 50/60 Hz, 100 Watts, 0.83 Amps
2050-1CEQ:	240 VAC, 50/60 Hz, 100 Watts, 0.42 Amps
2050-1CESUPQ:	240 VAC, 50/60 Hz, 100 Watts, 0.42 Amps
2050-1CEMARQ:	240 VAC, 50/60 Hz, 100 Watts, 0.42 Amps
2050-1CEIDQ:	240 VAC, 50/60 Hz, 100 Watts, 0.42 Amps
2052Q:	120 VAC, 50/60 Hz, 150 Watts, 1.25 Amps
2052IDQ:	120 VAC, 50/60 Hz, 150 Watts, 1.25 Amps
2052-1CEQ:	240 VAC, 50/60 Hz, 150 Watts, 0.63 Amps
2052-1CEIDQ:	240 VAC, 50/60 Hz, 150 Watts, 0.63 Amps
2053Q:	120 VAC, 50/60 Hz, 200 Watts, 1.67 Amps
2053-1CEQ:	240 VAC, 50/60 Hz, 200 Watts, 0.83 Amps
2054Q:	120 VAC, 50/60 Hz, 300 Watts, 2.50 Amps
2054IDQ:	120 VAC, 50/60 Hz, 300 Watts, 2.50 Amps
2054-1CEQ:	240 VAC, 50/60 Hz, 300 Watts, 1.25 Amps
2054-1CEIDQ:	240 VAC, 50/60 Hz, 300 Watts, 1.25 Amps
2056Q:	120 VAC, 50/60 Hz, 400 Watts, 3.60 Amps
2056-1CEQ:	240 VAC, 50/60 Hz, 400 Watts, 1.67 Amps

Unpacking and Installation

Shipping Carton

This should be inspected upon delivery. When received, carefully examine for any shipping damage before unpacking. If damage is discovered, the delivering carrier should both specify and sign for the damage on your copy of the delivery receipt.

Open the carton carefully making certain that all parts are accounted for before packaging materials are discarded. After unpacking, if damage is found, promptly report it to the carrier and request a damage inspection promptly.

IMPORTANT: Failure to request an inspection of damage within a few days after receipt of shipment absolves the carrier from any liability for damage: you must call for a damage inspection promptly.

Location

Find a convenient, draft-free location near a grounded electrical outlet meeting the power requirements listed on the unit nameplate.

The heater should be level, though operation does not depend on this, and away from edges of the laboratory bench where there is danger of accidental burns or knocking the unit over. There should be a space between the unit and any walls of 2 inches or more.

Heating Blocks

Select an aluminum-heating block designed for the test tubes or other vessels to be heated. Insert block(s) in the well on top of the unit.

Electrical Requirements

- Use a properly grounded electrical outlet of correct voltage and current handling capacity.
- Do not remove or modify grounded power plug. Use only properly grounded outlets to avoid shock hazard.



Note

Blocks have a threaded hole that serves a dual function: a) The furnished block puller is threaded into the hole to facilitate removal of blocks from the unit. b) The threaded hole can also be used for inserting a thermometer for temperature calibration procedure.



Note

Leave unit disconnected when not in use.

Operation

**Warning**

Do not use in the presence of flammable or combustible materials or explosive gases. Do not use in the presence of pressurized or sealed containers—fire or explosion may result, causing death or severe injury.

**Warning**

Do not heat any substance above a temperature which will cause it to emit toxic fumes—death or severe injury may result.

**Warning**

Do not add water to the heating block or the empty well of a unit connected to electrical power. This can result in an explosion with possibility of serious injury or death.

**Caution**

Turn the power switch to off and unplug unit when not in use. This extends the life of the heating element and eliminates the possibility of accidental burns.

Power Up

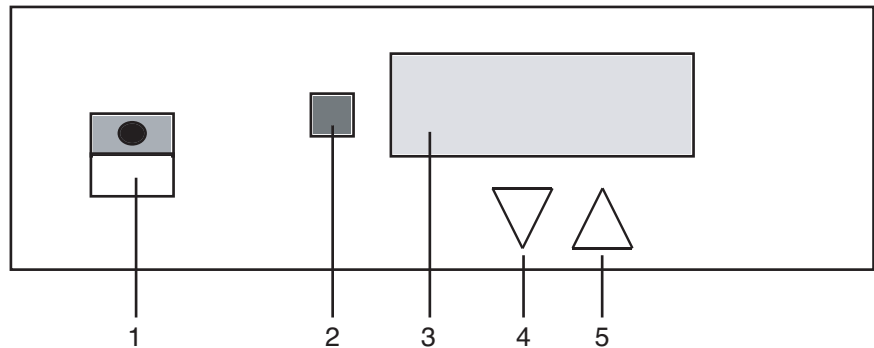
- Insert the plug into an outlet whose characteristics match those on the unit's data plate. Place the power switch in the ON position.

Positioning Block For Temperature Sensor Probe, 2000 Series

- Note the temperature sensor probe in the heater well—the bottom of the block has a hole that accommodates the sensor probe. Carefully insert the sensor probe into this hole in the bottom of the block.
- Insert test tubes or other vessels into the heater block once the temperature has stabilized at the desired point.

Note: To achieve the desired reaction or result, do not place materials in block until required temperature has been reached.

Control Panel: 2000 Series



1. **POWER SWITCH:** Turns the unit off and on.
2. **HEAT STATUS LAMP:** This lamp is lit when the heater is operating.
3. **TEMPERATURE DISPLAY:** Shows current operating temperature and set point when selected.
4. **UP ARROW KEY:** Press to increase set point.
5. **DOWN ARROW KEY:** Press to decrease set point.

Setting the Temperature: 2000 Series

- Turn the power switch ON. The display will flash the current revision number for about 5 seconds, then the temperature will be displayed.
- Press and release either the UP or DOWN ARROW KEY once—the display will flash the existing set point temperature already established.
- To change a temperature set point, press the appropriate UP or DOWN key to raise or lower the temperature to a desired value and release. When the displayed temperature stops flashing—showing the actual temperature—the new set point is established.

**Note**

If the up or down arrow keys are not touched for 5 seconds the display will default to the actual temperature.

**Note**

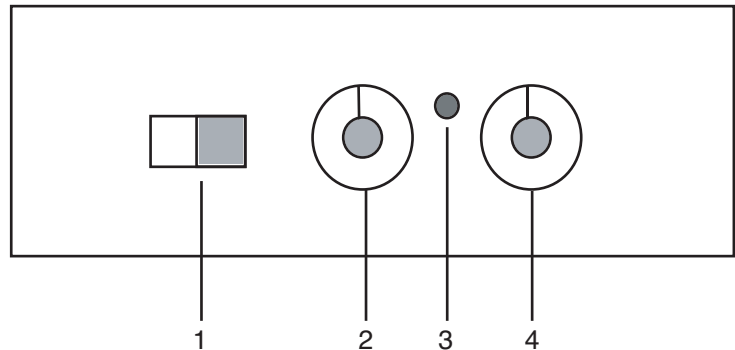
It is best to turn the power switch to off and unplug unit when not in use—this extends the life of the heating element and eliminates the possibility of accidental burns.

Temperature Calibration: 2000 Series

If the displayed temperature does not match the actual block temperature, calibrate the readout as follows:

- Press and hold both the UP and DOWN ARROW KEYS together at the same time until the display begins to flash.
- Press the appropriate UP or DOWN ARROW KEY to adjust the displayed temperature to match the actual temperature. Upon releasing the ARROW KEY, the display will stop flashing and the displayed temperature will match the thermometer temperature. Calibration data is automatically entered and stored.

Control Panel: 2050 Series



1. **POWER SWITCH:** Turns the unit off and on.
2. **LOW RANGE THERMOSTAT CONTROL:** Controls lower temperatures.
3. **HEAT STATUS LAMP:** This lamp is lit when the heater is operating.
4. **HIGH RANGE THERMOSTAT CONTROL:** Controls higher temperatures.

Setting the Temperature: 2050 Series

- Insert the plug into its outlet. Move the power switch from the OFF position either to the HIGH or the LOW position to select the desired heating range. The two ranges overlap: Low Range, from slightly above ambient to 60°C; High Range, from 50°C to 130°C. For an operating temperature between 50°C and 60°C, either thermostat can be used. The heater lamp will be lit when power is being supplied to the heating element.
- Set the thermostat control knob corresponding to the range selected. Dials are marked 0 to 10 for reference purposes only; the markings are not calibrated in degrees. As a rule of thumb, 0 on the low thermostat control indicates a

temperature slightly above ambient and 10 heats to about 60°C. On the high thermostat control, 0 represents a temperature of about 50°C and a setting of 10 produces about 130°C. A midpoint setting on the high thermostat control should produce a temperature of approximately 90°C.

- Check the thermometer. When the temperature has stabilized, make adjustments to raise or lower temperature as required. A flickering heater lamp indicates that the thermostat is maintaining a given temperature—it DOES NOT signal a malfunction.

Maintenance



Note

Make no attempt to service or repair a Thermo Scientific product under warranty before consulting your Thermo Scientific dealer. After the warranty period, such consultation is still advised, especially when the repair may be technically sophisticated or difficult. If assistance is needed beyond what the distributor can provide, please call Customer Service at 800-553-0039. No merchandise should be returned directly to the factory without obtaining a Return Materials Authorization (RMA) number from Customer Service.



Warning

Disconnect unit from power source before performing any service or repair.



Note

Users are solely responsible for the decontamination of their unit(s) if hazardous material is spilled on or into the unit(s).



Note

Before using any cleaning or decontamination method except those recommended by the manufacturer, users should check with customer service that the proposed method would not damage the equipment.

Since the dri-baths have no mechanical moving parts, they require very little maintenance beyond normal cleanup with a mild detergent and water.

- Turn the power switch to OFF and unplug the unit from its power source before attempting to clean up any spills or to perform maintenance, repairs or service. Allow unit to cool to near room temperature before removing the heating block(s).
- Remove the thermometer, where applicable, test tubes and/or other vessels from the heating block and wash them thoroughly or wipe them clean. Take the heating block from the well and wash in hot soapy water with a soft cloth. Holes in a block can be cleaned using a test tube brush. Rinse and wipe dry. Heating blocks can be submerged in hot water, but **UNDER NO CIRCUMSTANCES SHOULD THE CABINET BE IMMERSSED**. Wipe the exterior and well with a wet cloth, taking care not to get water in cabinet.
- Allow the cabinet to dry completely before installing blocks and reconnecting the power source.

Troubleshooting

The following is intended as a guide to help in servicing this unit, if problems should occur.



Note

Before attempting any repair, disconnect power cord from outlet. Depending upon controller options, not all of the following will relate to all models.

Symptom

Possible Cause of Problem

The heater status lamp will not light and the unit will not heat:

- Setpoint adjusted below actual temperature – readjust.
- Thermal cutoff is blown – replace.
- The unit is not plugged in, or it is not plugged into a proper, working outlet – check the outlet.
- The power switch is OFF – turn it ON.

Temperature is too high or too low and will not change when temperature setpoint is changed:

- Faulty temperature-control printed circuit board or sensor – check sensor resistance (110 | @ 25°C) and replace faulty component.
- Heater has failed – replace.
- Faulty thermostat – replace.
- Thermal cutoff is open – replace.

The heater status lamp flashes intermittently:

NOT A PROBLEM – the unit is turning itself on and off to maintain a constant temperature.

The heater status lamp is flicking but the unit will not heat to the desired level:

- The temperature control adjustment is set too low – turn clockwise.
- Faulty thermostat – replace.

The heater status lamp is lit, but not flickering; the unit is not hot:

- Allow more time to heat up.
- Faulty heater – replace.
- Thermal cutoff blown – replace.

Replacement Parts

DESCRIPTION	PART NUMBER		
Block Puller:	566-336-00	200 Watt, 240V, Model	
Cordset:		2002-1CEQ:	340-325-00
(120V)	CRX72	300 Watt, 120V, Model 2003Q:	340-331-00
(240V)	CRX70	300 Watt, 240V, Model	
Power Switch, 120 VAC:	440-359-00	2003-1CEQ:	340-326-00
Power Switch, 240 VAC:	440-292-00	400 Watt, 120V, Model 2004Q:	340-322-00
Rubber Feet (4):	790-214-00	400 Watt, 240V, Model	
RTD Temperature Probe,		2004-1CEQ:	340-322-01
2000 Series:	410-632-00	100 Watt, 120V, Model 2050Q:	340-330-00
Power Entry Module	CEX239	240V, Model 2050-1CEQ:	340-328-00
		150 Watt, 120V, Model 2052Q:	340-329-00
Fuses:		240V, Model 2052-1CEQ:	340-327-00
2000Q (2) 1.6aT	FZX47	200 Watt, 120V, Model 2053Q:	340-324-00
2002Q, 2003Q, (2) 4aT	FZX53	240V, Model 2053-1CEQ:	340-325-00
2004Q (2) 6.3aT	FZX61	300 Watt, 120V, Model 2054Q:	340-331-00
(2) 2aT, Models 2000-1CEQ,		240V, Model 2054-1CEQ:	340-326-00
20001Q, 2001-1CEQ, 2002-1CEQ,		400 Watt, 120V, Model 2056Q:	340-174-00
2003-1CEQ, 2004-1CEQ,		240V, Model 2056-1CEQ:	340-179-00
2050-1CEQ, 2052-1CEQ,		Electrical Reference Drawing,	
2053-1CEQ, 2054-1CEQ,		Model 2050:	228-692-00
2056-1CEQ:	5120-0025		
Fuse Drawer	FZX56	Wiring Schematics:	
		Models 2000Q~2004Q:	LT1873X2
Temperature Controller:		Models 2000-1CEQ~2004-1CEQ:	LT1873X3
2000Q, 2000-1CEQ, 2001Q,		2050Q, 2052Q, 2053Q, 2054Q,	
2001-1CEQ	PCX123	2056Q:	228-973-00
2002Q, 2002-1CEQ:	PCX124	2050-1CEQ, 2052-1CEQ,	
2003Q, 2003-1CEQ, 2004Q,		2053-1CEQ, 2054-1CEQ,	
2004-1CEQ:	PCX125	2056-1CEQ:	LT1960X1
Thermal Cutoff, 216°C:	330-236-00		
Power Switch (3-position),			
Model 2050:	440-373-00		
Status Lamp Base:	360-233-01		
Amber, Status Lamp Lens:	360-235-00		
Thermostat, High Range,			
Model 2050:	920-418-00		
Thermostat, Low Range,			
Model 2050:	920-417-00		
Thermostat Knob, order as set,			
Model 2050:	560-242-00		
Heaters:			
100 Watt, 120V, Model 2000Q:	340-330-00		
100 Watt, 240V, Model			
2000-1CEQ:	340-328-00		
150 Watt, 120V, Model			
2001Q:	340-329-00		
150 Watt, 240V Model			
2001-1CEQ:	340-327-00		
200 Watt, 120V, Model			
2002Q:	340-324-00		

Accessories

MODULAR BLOCK ACCESSORIES:

HEATING BLOCKS	NO. OF HOLES	CATALOG NO.
For 0.2 and 0.5ml tubes	19(0.2), 30(0.5)	2058Q
For 0.2ml tubes	79	2059Q
For 1 titer plate, size of 2 blocks*	0	2064Q
For 1 titer plate, size of 2 blocks	0	2065Q
For 10mm cuvettes	12	2066Q
For 8mm microcentrifuge tubes	30	2068Q
For 10mm microcentrifuge tubes	20	2069Q
For 6mm test tubes	30	2070Q
For 10mm test tubes	24	2071Q
For 12mm & 13mm test tubes	22	2072Q
For 15mm & 16mm test tubes	12	2073Q
For 20mm test tubes	8	2074Q
For 25mm test tubes	6	2075Q
3 Holes for 25mm test tubes,		
5 Holes for 12mm & 13mm test tubes		
6 Holes for 6mm test tubes	14	2076Q
For use as a hotplate	0	2078Q
For 17 and 18mm test tubes	12	2081Q
For 1 titer plate or 96 0.2 ml tubes and 6 0.5 ml tubes		2083Q
For 15 ml conical bottom centrifuge tubes	12	BKX40LLQ
For 50 ml conical bottom centrifuge tubes	4	BKX43LLQ
Stainless steel cover for Models 2064 and 2065		2085Q

*Direct Contact

Ordering Procedures

Please refer to the Specification Plate for the complete model number, serial number, and series number when requesting service, replacement parts or in any correspondence concerning this unit.

All parts listed herein may be ordered from the **Thermo Scientific** dealer from whom you purchased this unit or can be obtained promptly from the factory. When service or replacement parts are needed we ask that you check first with your dealer. If the dealer cannot handle your request, then contact our Customer Service Department at 1-866-9-thermo.

Prior to returning any materials, please contact our Customer Service Department for a "Return Materials Authorization" number (RMA). Material returned without an RMA number will be refused.

Two Year Limited Warranty

This Thermo Scientific product is warranted to be free of defects in materials and workmanship for two (2) years from the first to occur of (i) the date the product is sold by the manufacturer or (ii) the date the product is purchased by the original retail customer (the "Commencement Date"). Except as expressly stated above, the MANUFACTURER MAKES NO OTHER WARRANTY, EXPRESSED OR IMPLIED, WITH RESPECT TO THE PRODUCTS AND EXPRESSLY DISCLAIMS ANY AND ALL WARRANTIES, INCLUDING BUT NOT LIMITED TO, WARRANTIES OF DESIGN, MERCHANT ABILITY AND FITNESS FOR A PARTICULAR PURPOSE.

An authorized representative of the manufacturer must perform all warranty inspections. In the event of a defect covered by the warranty, we shall, as our sole obligation and exclusive remedy, provide free replacement parts to remedy the defective product. In addition, for products sold within the continental United States or Canada, the manufacturer shall provide free labor to repair the products with the replacement parts, but only for a period of ninety (90) days from the Commencement Date.

The warranty provided hereunder shall be null and void and without further force or effect if there is any (i) repair made to the product by a party other than the manufacturer or its duly authorized service representative, (ii) misuse (including use inconsistent with written operating instructions for the product), mishandling, contamination, overheating, modification or alteration of the product by any customer or third party or (iii) use of replacement parts that are obtained from a party who is not an authorized dealer of Thermo Scientific products.

Heating elements, because of their susceptibility to overheating and contamination, must be returned to the factory and if, upon inspection, it is concluded that failure is due to factors other than excessive high temperature or contamination, the manufacturer will provide warranty replacement. As a condition to the return of any product, or any constituent part thereof, to the factory, it shall be sent prepaid and a prior written authorization from the manufacturer assigning a Return Materials Number to the product or part shall be obtained.

IN NO EVENT SHALL THE MANUFACTURER BE LIABLE TO ANY PARTY FOR ANY DIRECT, INDIRECT, SPECIAL, INCIDENTAL, OR CONSEQUENTIAL DAMAGES, OR FOR ANY DAMAGES RESULTING FROM LOSS OF USE OR PROFITS, ANTICIPATED OR OTHERWISE, ARISING OUT OF OR IN CONNECTION WITH THE SALE, USE OR PERFORMANCE OF ANY PRODUCTS, WHETHER SUCH CLAIM IS BASED ON CONTRACT, TORT (INCLUDING NEGLIGENCE), ANY THEORY OF STRICT LIABILITY OR REGULATORY ACTION.

For the name of the authorized Thermo Scientific product dealer nearest you or any additional information, contact us:

308 Ridgefield Court, Asheville, NC, 28806 USA

Phone: 1-866-984-3766

Fax: 1-828-665-4071

Web: www.thermo.com