## Thermo Scientific System III Heat Exchanger

Thermo Scientific Manual P/N U00678 Rev. 09/28/2015

## Installation Operation Basic Maintenance

CE

Visit our Web site at:

http://www.thermoscientific.com/tc Product Service Information, Applications Notes, MSDS Forms, e-mail.

Voice Info: (800) 258-0830



### Thermo Scientific System III Heat Exchanger

Thermo Scientific Manual P/N U00678 Rev. 09/28/2015

## Installation Operation Basic Maintenance

CE

Visit our Web site at:

http://www.thermoscientific.com/tc Product Service Information, Applications Notes, MSDS Forms, e-mail.

Voice Info: (800) 258-0830

#### **Thermo Fisher Scientific**

25 Nimble Hill Road Newington, NH 03801 Tel : (800) 258-0830 or (603) 436-9444 Fax : (603) 436-8411 www.thermoscientific.com/tc

#### Sales, Service, and Customer Support

25 Nimble Hill Road Newington, NH 03801 Tel: (800) 258-0830 Sales: 8:00 am to 5:00 pm Service and Support: 8:00 am to 6:00 pm Monday through Friday (Eastern Time) Fax: (603) 436-8411 service.tc.us@thermofisher.com

Dieselstrasse 4 D-76227 Karlsruhe, Germany Tel : +49 (0) 721 4094 444 Fax : +49 (0) 721 4094 300 info.tc.de@thermofisher.com

Building 6, No. 27 Xin Jinqiao Rd., Shanghai 201206 Tel : +86(21) 68654588 Fax : +86(21) 64457830 info.china@thermofisher.com

#### Statement of Copyright

Copyright © 2015 Thermo Fisher Scientific. All rights reserved. This manual is copyrighted by Thermo Fisher Scientific. Users are forbidden to reproduce, republish, redistribute, or resell any materials from this manual in either machinereadable form or any other form.

## System III Liquid to Liquid Heat Exchanger Instruction and Operation Manual

### **Table of Contents**

Preface		
	Compliance	3
	WEEE/RoHS	3
	After-sale Support	3
	Unpacking	4
	Warranty	4
	Feedback	4
Section I Safety		
	Warnings	5
Section II General Information		
	Description	7
	Specifications	7
Section III Installation		
	Site	
	Facility Water Requirements	
	Electrical Requirements	
	Voltage Selection	
	Plumbing Requirements	
	External Pressure Regulator (Optional)	
	Fluids	
	Filling Requirements	
	Automatic Refill Device (Optional)	
	Flow Control	
		12
Section IV Operation		
	Start Up	
	Temperature Controller	
	Setup Loop	15

### Section V Maintenance& Service

Service Contracts	17
Pump Strainer	17
Facility Water Strainer	18
Algae	18
Configuration	18
Voltage Selection	18
Pressure Relief Valve	19
Fault Interlock Contact	19
Fuses	19
Phase Rotation	19
Pump Motor Overload Protector	20
Displaying Software Version	20

### Section VI Troubleshooting

Error Codes	21
Poor Temperature Stability	22
Service Assistance	22
Parts List	22

### Section VII Diagrams

TU Pump Flow Diagram	23
CP Pump Flow Diagram	24
Front View	
Rear View	25
Large Volume Exchangers	26

### Appendix

### Water Quality and Standards

### WARRANTY

### Preface

#### Compliance

Refer to the Declaration of Conformity that shipped with the exchanger.

#### WEEE/RoHS

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. 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 Thermo Fisher Scientific's compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at: **www.thermofisher.com/WEEERoHS** 

### **After-sale Support**

Thermo Fisher Scientific is committed to customer service during and after the sale. If you have questions concerning the operation of your exchanger, contact our Sales Department. If your exchanger fails to operate properly, or if you have questions concerning spare parts, contact our Customer Service Department. Before calling, please obtain the following information:

- BOM number
- Serial number
- Software version

The BOM and serial number are on a label on the rear of the exchanger. To display the software version see page 20.

#### Unpacking

Retain all cartons and packing material until the exchanger has been operated and found to be in good condition. If the exchanger shows external or internal damage, contact the transportation company and file a damage claim. Under ICC regulations, this is your responsibility.

#### Out of Box Failure

An out of box failure is defined as any product that fails to operate in conformance with the seller's published specifications at initial power up. The exchanger must be installed in accordance with manufacturer's recommended operating conditions within 30 days of shipment from the seller.

Any product meeting the definition of an out of box failure must be packed and shipped back in the original packaging to Thermo Fisher Scientific for replacement with a new exchanger. The seller pays all shipping costs. The customer must obtain a Return Material Authorization (RMA) from Thermo Fisher prior to shipping.

#### Warranty

Exchangers have a warranty against defective parts and workmanship for one full year from date of shipment. See back page for more details.

#### Feedback

We appreciate any feedback you can give us on this manual. Please e-mail us at

tcmanuals@thermofisher.com.

Be sure to include the manual part number and the revision date listed on the front cover.

### **Section I Safety**

#### Warnings

Make sure you read and understand all instructions and safety precautions listed in this manual before installing or operating your exchanger. If you have any questions concerning the operation of your exchanger or the information in this manual, contact our Sales Department.

Performance of installation, operation, or maintenance procedures other than those described in this manual may result in a hazardous situation and may void the manufacturer's warranty.

Observe all warning labels.

Never remove warning labels.

Never operate damaged or leaking equipment.

Always turn off the exchanger and disconnect the line cord from the power source before performing any service or maintenance procedures, or before moving the exchanger.

Always empty the reservoir before moving the exchanger.

Never operate equipment with damaged line cords.

Never operate without fluid.

Refer service and repairs to a qualified technician.

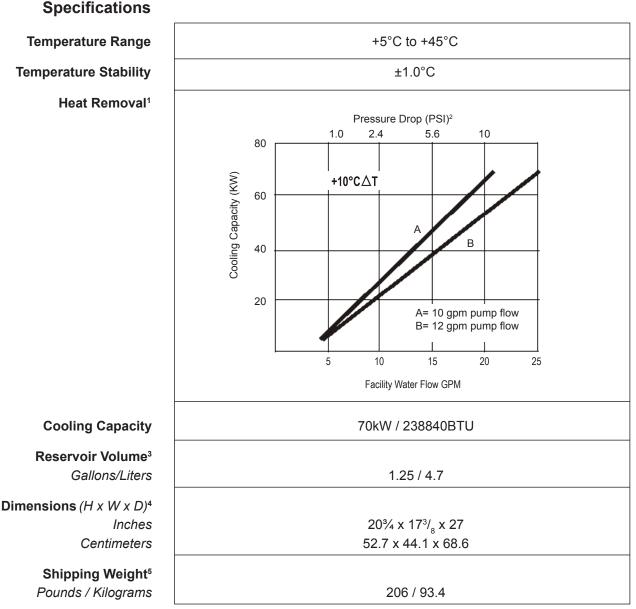


In addition to the safety warnings listed above, warnings are posted throughout the manual. These warnings are designated by an exclamation mark inside an equilateral triangle with text highlighted in bold print. Read and follow these important instructions. Failure to observe these instructions can result in permanent damage to the exchanger, significant property damage, or personal injury or death.

### **Section II General Information**

#### Description

The Thermo Scientific System III Liquid to Liquid Heat Exchanger is designed to remove heat from water-cooled instruments. The heat exchanger consists of a recirculation pump, PVC reservoir, and a microprocessor controller.



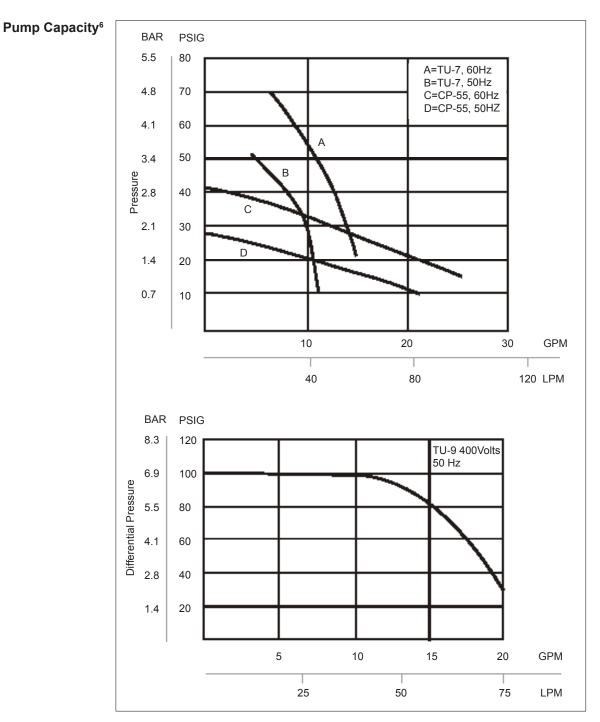
1. Specifications obtained using water as the recirculating fluid and as a coolant on the secondary loop, +25°C ambient condition, at nominal operating voltage. Other fluids, process temperatures, ambient temperatures, altitude, or operating voltages will affect performance. Specifications subject to change. Heat load removal is based on a 10°C  $\Delta$  between the temperature of the cooling water supply (house water) and the cooling fluid as it leaves the System. 2. Pressure drop obtained with the modulating valve fully open. Pressure drop indicates the minimum pressure differential pressure between the facility water inlet and outlet to achieve the

corresponding facility water flow rate (Pressure<sub>inlet</sub> - Pressure<sub>outlet</sub> = Pressure<sub>drop</sub>).

3. Larger volume reservoir exchangers are available.

4. Exchangers with CP 55 pumps. Exchangers with TU 7 and TU 9 pumps measure 33<sup>3</sup>/<sub>4</sub> x 23 x 27 (85.7 x 58.4 x 68.6). For complete dimensions see pages 25-26.

5. Larger volume exchangers weigh approximately 289 pounds (131 kilograms).



6. TU pumps require 3Ø power. TU 7 and TU 9 requires large case.

### Section III Installation

#### Site

The exchanger should be located in a laboratory or clean industrial environment with easy access to a facility cooling water supply and a drain. Never place the exchanger in a location where excessive heat, moisture, or corrosive materials are present.

Refer to the pump label on the rear of the exchanger to identify the specific type of pump. Exchangers with a TU-7 pump are equipped with a pump motor fan. The fan is used to cool the pump motor and prevent the motor from overheating. Air is drawn through the front of the exchanger and is discharged through the rear. A minimum clearance of 6 inches (0.15 meters) at the front and rear of the exchanger is necessary for ventilation.

### Facility Water Requirements



Limit the facility water inlet pressure to less than 150 psi (10,2 Bar) and limit the facility water inlet pressure to outlet pressure differential across the System III to less than 35 psid (2,4 Bar).

Refer to the Cooling Capacity chart in Section II, Specifications. The flow rate of the cooling water supply must meet or exceed these requirements for the exchanger to operate at its full rated capacity. If the cooling water does not meet these standards, the cooling capacity will be derated. The chart is based on a difference between the temperature of the cooling water supply (House Water) and the cooling fluid flowing from the System III to the instrument being cooled.

As the heat load increases, the required flow rate of the cooling water supply increases. For example, on a System III with a 6 gpm pump flow, if the heat load is 12 kilowatts, approximately 3 gpm of cooling water flow is required to remove the heat. However, if the heat load is increased to 36 kilowatts, about 8 gpm of cooling water flow is required.

The flow meter on the front of the exchanger does not measure the flow rate of the cooling water supply. The flow meter measures the flow rate of the cooling fluid going to the instrument being cooled.

### Electrical Requirements



The exchanger construction provides protection against the risk of electric shock by grounding appropriate metal parts. The protection will not function unless the power cord is connected to a properly grounded outlet. It is the user's responsibility to assure a proper ground connection is provided.

Refer to the serial number label on the rear of the exchanger for the specific electrical requirements.

Ensure the voltage of the power source meets the specified voltage, ±10%.

Pump	Voltage	Frequency	Phase	Amperage	Line Cord Plug
CP55	230	60	1	9	L6-15
TU7	230	60	3	8.2	L15-20
TU9	230	60	3	13.8	L15-20
TU7	380	50	3 (5W)	4	None

# Voltage Selection (60 Hertz Exchangers)

If the exchanger is to be operated from a 220 to 240V source, a voltage range selector switch inside the exchanger must be reset. See Maintenance and Service section for instructions on changing the voltage selector.

NOTE Exchangers rated 380-415 volt do not have a voltage selection switch.

### Plumbing Requirements

Before installing the exchanger to an instrument that previously used tap water as a cooling fluid, flush the instrument several times to remove any rust or scale that has built up. The manufacturer of the instrument should be able to recommend a cleaning fluid for their equipment.

The plumbing connections are located on the rear of the exchanger and are labelled FACILITY WATER and RECIRCULATING CLEAN FLUID.

The connections for exchangers with CP pumps are all are all 1 inch FPT. The facility water connections for exchangers with TU pumps are 1 inch, the process fluid connections are 1 1/2 inch.

A basket strainer is supplied with the exchanger to protect the heat exchanger from becoming clogged by dirty cooling water. Install this strainer on the FACILITY WATER inlet. A clogged strainer can adversely affect cooling capacity. See Section V, Facility Water Strainer for cleaning instructions.

Connect the FACILITY WATER connections to the cooling water supply and the drain. Connect the RECIRCULATING CLEAN FLUID connections to the instrument being cooled.

Flexible tubing, if used, should be of heavy wall or reinforced construction. All tubing should be rated to withstand 80 psi at +40°C. Make sure all tubing connections are securely clamped. Avoid running tubing near radiators, hot water pipes, etc. If substantial lengths of tubing are necessary, insulation may be required to prevent the loss of cooling capacity.

Tubing and insulation are available from Thermo Fisher. Contact our Sales Department for more information (see Preface, After-sale Support).

It is important to keep the distance between the exchanger and the instrument being cooled as short as possible, and to use the largest diameter tubing practical. Tubing should be straight and without bends. If reductions must be made, they should be made at the inlet and outlet of the instrument being cooled, not at the exchanger.

### External Pressure Regulator (Optional)

An optional External Pressure Reducer (EPR) limits the maximum pressure to your application. We recommend an EPR when circulating to applications that are sensitive to higher pressures.

### Fluids



Never use flammable or corrosive fluids. Do not use automotive antifreeze. Commercial anti-freeze contains silicates that can damage the pump seals. Use of automotive anti-freeze will void the manufacturer's warranty.

We recommend distilled/deionized water with a 0.05 - 0.1 MOhmcm reading.



Highly distilled/deionized water, above the 3 MOhmcm region, may become aggressive and is not recommended for use with exchangers with wetted parts other than stainless steel. Distilled/deionized water in the 15 MOhmcm region is definitely aggressive and should not be used. Exchangers operating in these regions should be closely monitored. See Water Quality Standards and Recommendations in the Appendix.

If you do not have access to distilled/deionized water we recommend using filtered tap water. Thermo Fisher cannot recommend any custom fluids, these fluids are too dependent on your particular application.

Below +8°C a non-freezing solution is required. The selected cooling fluid must have a viscosity of 50 centistokes or less. A 50/50 mixture, by volume, of distilled/deionized water and laboratory grade ethylene glycol is suggested.

Filling Requirements	
Requirements	If substantial lengths of cooling lines are required, they should be pre-filled with cooling fluid before connecting them to the exchanger.
	Remove the reservoir cover. Fill the reservoir with clean cooling fluid to within one inch of the top of the reservoir. Have extra cooling fluid on hand and follow the steps described in Section IV, Start Up.
Automatic Refill Device (Optional)	
	The automatic refill device maintains the correct level of cooling fluid in the reservoir. The device consists of a float switch in the reservoir and a sole- noid valve at the rear of the exchanger. If the cooling fluid level falls, the float switch will drop, opening the solenoid valve and allowing make-up fluid to fill the reservoir. Once the cooling fluid reaches the proper level, the float switch will rise and the solenoid valve will close.
	<b>NOTE:</b> For the solenoid valve to close properly, the minimum supply flow rate should be 2.2 gallons per minute.
	Connect the $3/_{8}$ inch OD stainless steel barbed fitting on the solenoid value to the make-up fluid source using $5/_{16}$ or $3/_{8}$ inch ID flexible tubing.
	Tubing is available from Thermo Fisher. Contact our Sales Department for more information (see Preface, After-sale Support).
Flow Control	The RECIRCULATING FLOW CONTROL handle is connected to a three-way valve that controls the flow of the cooling fluid to the instrument being cooled. The handle is located on the front of the exchanger.
	When the handle is in the "+" position, the valve is open and all available cooling fluid is supplied to the instrument being cooled. When the handle is in the " <b>0</b> " position, the valve is closed and no cooling fluid is supplied to the instrument being cooled. When the handle is between these two positions, the flow rate of the cooling fluid is between full flow and no flow. Use the flow meter to adjust the desired flow rate.
	The gauge next to the flow control handle indicates the operating pressure.
Auto Restart	Exchangers are equipped with an auto restart feature. If power is lost, the

Exchangers are equipped with an auto restart feature. If power is lost, the exchanger will automatically restart when power is restored. This feature is enabled/disabled using the controller's Setup Loop, see page 15.

### **Section IV Operation**

### Start Up

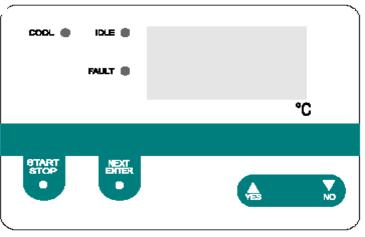
Before starting, double check all electrical and plumbing connections and make sure the circulating system (the System III, your application, and the tubing that connects them) has been properly filled with cooling fluid.

Turn the RECIRCULATING FLOW CONTROL to "**0**". Press START/STOP. The pump starts and the controller displays the reservoir fluid temperature.

The low fluid level monitor in the reservoir prevents the exchanger from operating if the fluid level in the reservoir is below a safe level. By slightly, and/or intermittently opening the RECIRCULATING FLOW CONTROL (toward "+") and using extra cooling fluid to keep the reservoir topped off, the system can be filled without repeated tripping of the low fluid level monitor.

If the exchanger shuts down, top off the reservoir and restart it. When the system is full, the reservoir level will no longer drop when the RECIRCULAT-ING FLOW CONTROL valve is opened (toward "+").

A control valve, located in the FACILITY WATER inlet line, regulates the flow rate of the cooling water supply as it enters the exchanger. The valve regulates the flow rate based on the heat load. Flow through the exchanger stops automatically when the exchanger is shut off.



Temperature Controller

To turn the exchanger off, press START/STOP. The pump will stop.

The IDLE and COOL lights indicate the control valve's status. When the temperature control valve is wide open (for maximum cooling), the COOL light is on steady. When the control valve is closed, the IDLE light is on. As the control valve moves between these extremes, the two lights flash with varying on-time to indicate the approximate position of the control valve.

#### **Temperature Controller**

The microprocessor controller controls temperature using a PID (Proportional-Integral- Derivative) algorithm. It is designed with self-diagnostic features and easy to use operator interface.

#### NEXT ENTER

Use this key to accept and save changes.

### YES, 📥

This key is used to increase numerical values.

### NO, 🔻

This key is used to decrease numerical values.

When the controller is powered it displays the reservoir fluid temperature. Press the **NEXT ENTER** key to view the setpoint. The display flashes between **SP** and the actual setpoint number. If desired, use the **YES** and **NO** keys to change the setpoint. The display flashes as soon as either key is depressed. Once the desired setpoint is displayed, press **NEXT ENTER**.

When selecting an operating temperature, remember that the lowest achievable temperature is a function of the available flow rate, the temperature of the cooling water supply and the heat load.

The temperature control system actuates a control valve in the FACILITY WATER line. The control valve adjusts the flow of the cooling water supply to produce the desired operating temperature.

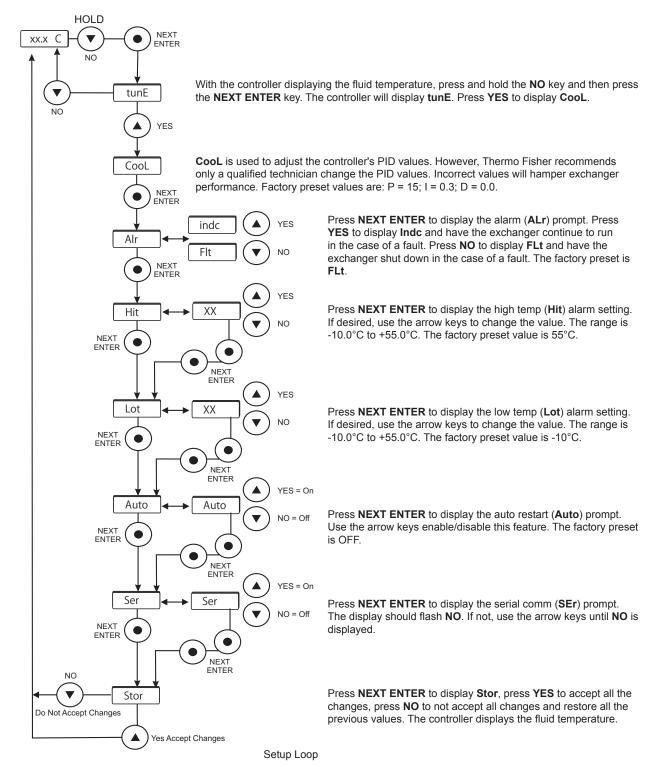
**NOTE:** The controller does not use the new value until the **NEXT ENTER** key is depressed and the display stops flashing. The controller will not allow you to enter a value above the maximum or below the minimum value, or any illegal value. If you try to enter an illegal value the display will revert to its original value when the last digit was entered.

If the **NEXT ENTER** key is not depressed within one minute, the controller will time out and the new value will not be accepted. The controller will revert to the previous value.

NOTE: Error codes are addressed in Section V, Maintenance and Service.

The controller is used to tune the COOL PID values; configure the exchanger to continue to run (Indc) or shut down (FLt) in the event of a fault; set the high (Hit) and low (Lot) temperature limits; and enable/disable auto restart (Auto). If a temperature limit is exceeded the controller will display an error code, see Section VI.

**NOTE:** Serial Communication (SEr) is not operational. Ensure the display indicates OFF.



### Section V Maintenance & Service



For personal safety and equipment reliability, the following procedure should only be performed by a qualified technician. Contact our Service Department for assistance (see Preface, After-sale Support).

#### Service Contracts

Thermo Fisher Scientific offers on-site Service Contracts that are designed to provide extended life and minimal down-time for your exchanger. For more information, contact our Service Department (see Preface, After-sale Support).

#### **Pump Strainer**

Exchangers with TU pumps have a wire mesh pump strainer is located at the bottom of the reservoir. If debris is drawn into the reservoir, the strainer will prevent the material from being sucked into the pump and damaging the pump vanes.

After initial installation, the strainer may become clogged with debris and scale within the first week. Therefore, the strainer must be cleaned after the first week of installation. After this first cleaning, the frequency of cleaning depends on the purity of the cooling water. We recommend a visual inspection of the reservoir be made monthly after the initial cleaning. After several months, the frequency of cleaning will be established.

If the strainer is visibly clogged, cleaning is required.

Disconnect the power cord from the power source and drain the reservoir before cleaning the strainer.

Cover the strainer with a plastic bag to help trap any particles which may become dislodged, and then remove the strainer by unscrewing it.

Clean the strainer by rinsing it with water.

Refer to Section III, Filling Requirements for instructions on replacing the cooling fluid.

### Facility Water Strainer

The facility water strainer is the user-installed basket strainer on the FACILITY WATER inlet. **NOTE:** The strainer is designed to be used only with water. Clean the strainer when it becomes visibly clogged or dirty.

Disconnect the power cord from the power source and turn off the facility cooling water.

Place a container under the strainer to collect any water that spills out of the basket when it is removed.

Unscrew the clear plastic basket. Remove the screen and rinse it with water. Replace the screen and the basket.

### Algae

To restrict the growth of algae in the reservoir, it is recommended that the reservoir cover be kept in place and that all circulation lines be opaque. This will eliminate the entrance of light which is required for the growth of most common algae.

### Configuration

The exchanger top is secured to the cabinet by four ball stud retainers; one at each per corner. Remove cabinet top by prying upward gently (cover pops off) in order to perform the following adjustments.

In some cases, the side access panels needs to be removed. The access panels are secured using screws installed through the bottom of the cabinet.

**NOTE**: Some models have a one piece cover rather than separate top and side panels. One piece covers are secured by five screws on each side of the case.

### Voltage Selection (60 Hertz Exchangers)

**NOTE:** Exchangers rated 380-415 volt do not have a voltage selection switch.

Remove the top cover from the cabinet, see Configuration.

The VOLTAGE SELECT toggle switch is located on the right side of the control box. Two ranges are available: 200-208V and 220-240V. Set the switch for the appropriate range.

Replace the top cover.

### Pressure Relief Valve

Exchangers with TU pumps have a pressure relief valve is located on the pump discharge line. The relief valve establishes the exchanger's maximum operating pressure. If the pressure of the fluid leaving the pump exceeds the valve setting, the relief valve will bypass the fluid within the exchanger to relieve to the pressure. The relief valve does not determine the actual operating pressure; the actual operating pressure is determined by the flow control valve setting and pressure drop through the instrument being cooled. If adjustment is necessary, consult our Service Department for assistance (see Preface, After-sale Support). Fault Interlock Contact A set of contacts is connected to a receptacle on the control panel. The contacts are rated 10A/240V. This is not a power inlet or outlet. The receptacle is isolated from the circuitry. Its ground pin is connected to the chassis. The contacts are closed during normal operation and open when the exchanger is turned off or when a fault is detected. **Fuses** 

Three-phase exchangers have fuses. Refer to the serial number label on the rear of the exchanger for the specific electrical requirements.

Remove the top cover, the right access panel and the control box cover. The fuses are located inside the control box.

### Phase Rotation

Three phase exchangers with three phase pump motors are equipped with a phase rotation interlock. If the phasing is wrong, the controller will display PHEr and the exchanger will not start.

Unplug the exchanger. Reverse any two power cord wires in the power cord plug.



#### Never remove the green ground wire.

Plug in the exchanger. The PHEr display should be off and the exchanger should start.

Pump Motor Overload Protector			
	Three phase exchangers with three phase pump motors have a pump motor overload protector.		
	The overload protector prevents the pump motor from exposure to excessively high current. If an overload fault occurs, due, for example, to a heavy work load, the controller will display <b>OL</b> and the exchanger will shut down. The overload protector will automatically reset after about two minutes. The exchanger must be manually restarted.		
Displaying Software Version			
	This procedure will display the software version number on the exchanger's display. In the event the exchanger is not operational, the software version (and the checksum) can also be read from the label which is on the micro-processor chip itself.		
	The following chart uses an example of software versi	on 000507.84A	
	begin at the reservoir temperature display	<b>20.3</b> °C	
	Press and hold NO for at least 10 seconds.	0507	
	displays software version digits to <b>left</b> of decimal. Note do not display.	e the two leading zeros	
	Press NEXT	84	
	displays software version digits to <b>right</b> of decimal.		
	Press NEXT	1	
	displays software version revision letter (as its equival cannot show letters. A=1, B=2, etc.)	ent number - display	
	Press NEXT	0000	
	displays checksum - this can be disregarded		
	Press NEXT	<b>20.3</b> °C	
	returns to reservoir temperature display		

### **Section VI Troubleshooting**

### Error Codes

The controller also has the capability to display error codes.

#### **Power up Errors**

Power up errors are displays until any controller key is pressed. If the error message persists the problem is likely a keypad or controller board failure.

Display	Indication
Er 00	ROM checksum
Er 01	RAM check
Er 02	Keypad failure
Er 03	NOVRAM checksum error

#### **Operating Errors**

Operating errors are displays once and then clear after three seconds or when any controller key is pressed. If error messages Er 04 - Er 14 persist, the problem is likely controller board failure. If Er 15 is displayed, ensure serial communications is turned off in the Setup Loop. If Er 16 is displayed the exchanger needs calibration, contact Thermo Scientific for assistance.

### Latching Errors

In order to restart the exchanger when a latching error is displayed, the controller's START/STOP button must be pressed after the condition is cleared.

Display	Indication
Lo t	Fluid temperature lower than low temp alarm setting.
Hi t	Fluid temperature higher than high temp alarm setting.
Er 22	Fluid temperature higher than the high overrange
	point. This is a fixed value, factory set at 55°C.
Er 25	RDT1 input shorted.
Er 26	RTD1 input open.
LLF	Low level fault switch open for at least three seconds.
PHEr	Improper phasing.
OL	Pump motor overload.

If any other code appears contact Thermo Fisher customer service, see Preface.

### Poor Temperature Stability

The System III is designed to maximize heat removal with minimal facility cooling water requirements. Under conditions of excessive cooling capacity, instability may result. The condition may be aggravated by : High temperature setpoint

Low facility water temperature

Small heat load

High facility water pressure

The situation can be remedied by limiting the available facility cooling water. Facility water pressure can be reduced with a pressure regulator before the System III. However, excessive restriction of facility water will reduce the System III's capacity.

### **Service Assistance**

If, after following these troubleshooting steps, your exchanger fails to operate properly, contact our Service Department for assistance (see Preface, Aftersale Support). Before calling please obtain the following information:

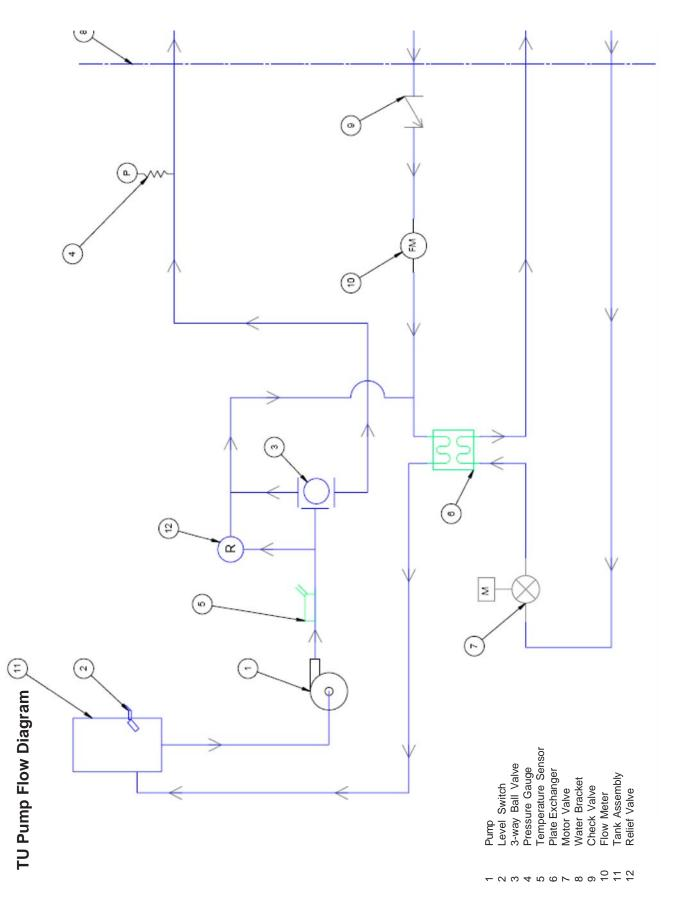
BOM number Serial number Voltage of exchanger Voltage of power source Temperature at which the problem occurs Temperature, pressure and flow rate of cooling water supply Software Version

### Parts List

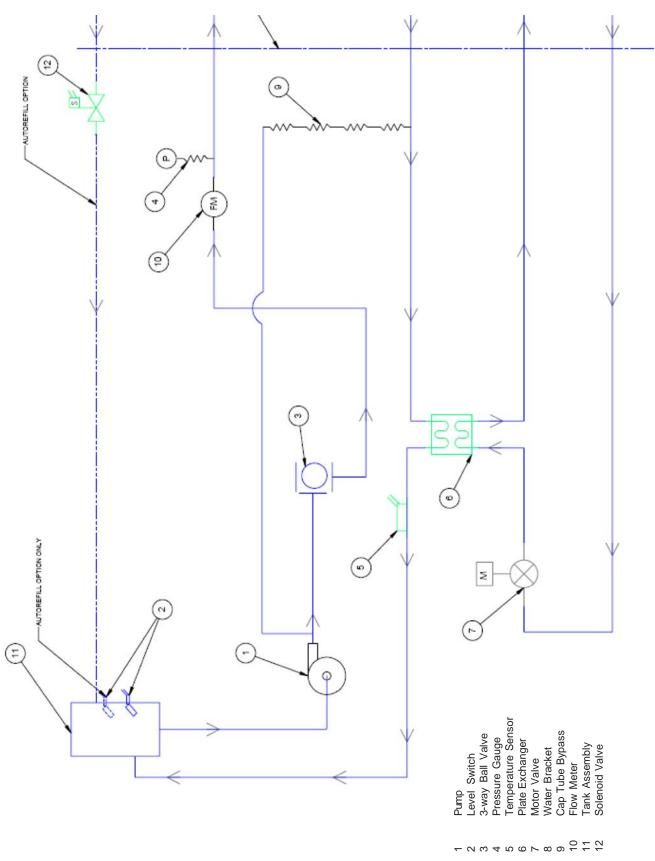
Our Service Department can provide you with a complete list of spare parts for your exchanger (see Preface, After-sale Support). Before calling, please obtain the following information:

BOM number

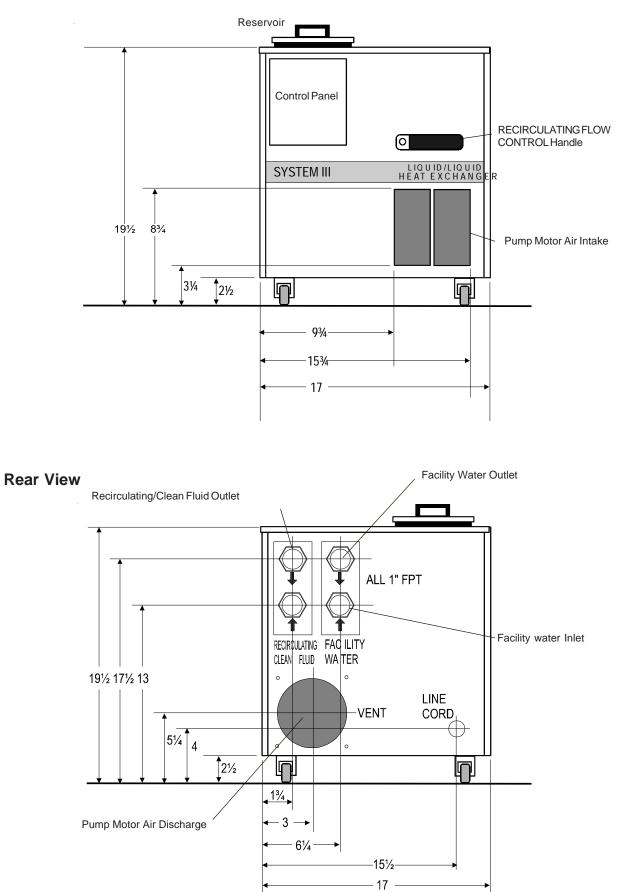
Serial number



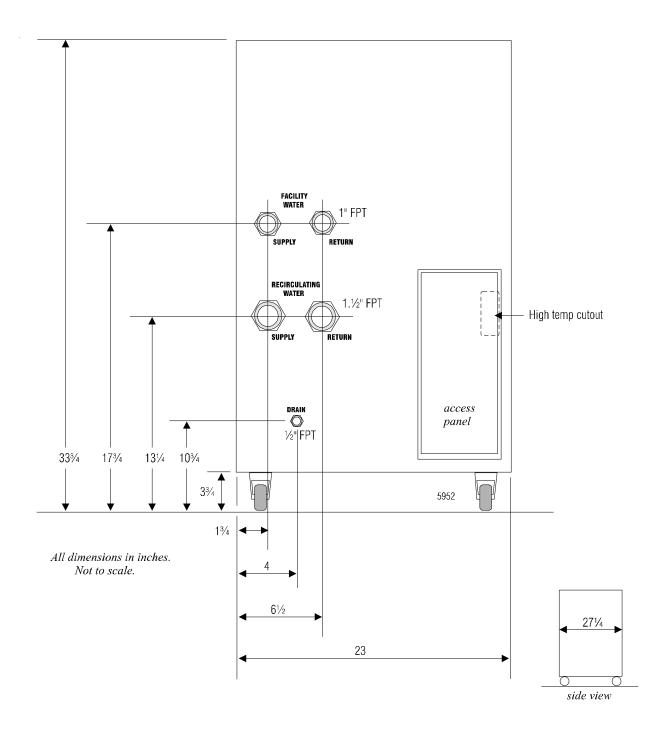




### **Front View**



### Large Volume Units



### Appendix

## Water Quality Standards and Recommendations

	Permissible (PPM)	Desirable (PPM)
Microbiological	S	
(algae, bacteria, fungi)	0	0
Inorganic Chemicals		
Calcium	<40	<0.6
Chloride	<250	<25
Copper	<1.3	<1.0
Iron	<0.3	<0.1
Lead	<0.015	0
Magnesium	<12	<0.1
Manganese	<0.05	< 0.03
Nitrates\Nitrites	<10 as N	0
Potassium	<20	<0.3
Silicate	<25	<1.0
Sodium	<20	<0.3
Sulfate	<250	<50
Hardness	<17	<0.05
Total Dissolved Solids	<50	<10
Other Parameters		
рН	6.5-8.5	7-8
Resistivity	0.01*	0.05-0.1*
* MOhmcm (Compensate	d to 25°C)	

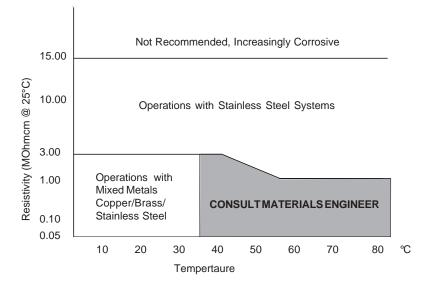
Unfavorably high total ionized solids (TIS) can accelerate the rate of galvanic corrosion. These contaminants can function as electrolytes which increase the potential for galvanic cell corrosion and lead to localized corrosion such as pitting. Eventually, the pitting could become so extensive that leaking will occur between the process water and facility water diminishing the System's heat transfer capability.

High water hardness (Calcium and Maganese) can also produce scaling. Scaling will inhibit heat transfer between the process and facility side by building up a deposit layer on metal surfaces. As an example, raw water in the United States averages 171 ppm (of NaCl). The recommended level for use in a water system is between 0.5 to 5.0 ppm (of NaCl).

Recommendation: Initially fill the tank with distilled/deionized water. Do not use untreated tap water as the total ionized solids level may be too high.

The desired level for long time usage is 1 to 3 MOhmcm (compensated to 25°C).

The above recommendation will reduce the electrolytic potential of the water and prevent or reduce the galvanic corrosion observed.



### WARRANTY

Thermo Fisher Scientific warrants for 12 months from date of shipment any Thermo Scientific product according to the following terms.

Any part of the product manufactured or supplied by Thermo Fisher Scientific and found in the reasonable judgment of Thermo Fisher to be defective in material or workmanship will be repaired at an authorized Thermo Fisher Repair Depot without charge for parts or labor. The product, including any defective part must be returned to an authorized Thermo Fisher Repair Depot within the warranty period. The expense of returning the product to the authorized Thermo Fisher Repair Depot for warranty service will be paid for by the buyer. Our responsibility in respect to warranty claims is limited to performing the required repairs or replacements, and no claim of breach of warranty shall be cause for cancellation or recision of the contract of sales of any product. With respect to products that qualify for field service repairs, Thermo Fisher Scientific's responsibility is limited to the component parts necessary for the repair and the labor that is required on site to perform the repair. Any travel labor or mileage charges are the financial responsibility of the buyer.

The buyer shall be responsible for any evaluation or warranty service call (including labor charges) if no defects are found with the Thermo Scientific product.

This warranty does not cover any product that has been subject to misuse, neglect, or accident. This warranty does not apply to any damage to the product that is the result of improper installation or maintenance, or to any product that has been operated or maintained in any way contrary to the operating or maintenance instructions specified in this Instruction and Operation Manual. This warranty does not cover any product that has been altered or modified so as to change its intended use.

In addition, this warranty does not extend to repairs made by the use of parts, accessories, or fluids which are either incompatible with the product or adversely affect its operation, performance, or durability.

Thermo Fisher Scientific reserves the right to change or improve the design of any product without assuming any obligation to modify any product previously manufactured.

THE FOREGOING EXPRESS WARRANTY IS IN LIEU OF ALL OTHER WARRANTIES, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO WARRANTIES OR MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE.

OUR OBLIGATION UNDER THIS WARRANTY IS STRICTLY AND EXCLUSIVELY LIMITED TO THE REPAIR OR REPLACEMENT OF DEFECTIVE COMPONENT PARTS AND Thermo Fisher Scientific DOES NOT ASSUME OR AUTHORIZE ANYONE TO ASSUME FOR IT ANY OTHER OBLIGATION.

Thermo Fisher ScientificASSUMES NO RESPONSIBILITY FOR INCIDENTAL, CONSEQUENTIAL, OR OTHER DAMAGES INCLUDING, BUT NOT LIMITED TO LOSS OR DAMAGE TO PROPERTY, LOSS OF PROFITS OR REVENUE, LOSS OF THE PRODUCT, LOSS OF TIME, OR INCONVENIENCE.

This warranty applies to products sold by Thermo Fisher Scientific. (Refer to the warranty for products sold by the affiliated marketing company of Thermo Fisher Scientific for any additional terms.) This warranty and all matters arising pursuant to it shall be governed by the law of the State of New Hampshire, United States. All legal actions brought in relation hereto shall be filed in the appropriate state or federal courts in New Hampshire, unless waived by Thermo Fisher Scientific.

Thermo Fisher Scientific 81 Wyman Street P.O. Box 9046 Waltham, Massachusetts 02454-9046 United States

www.thermofisher.com

