

Thermo Scientific System IV Heat Exchanger

Thermo Scientific Manual P/N U00682
Rev. 09/16/2015

Installation Operation Basic Maintenance



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 IV Heat Exchanger

Thermo Scientific Manual P/N U00682

Rev. 09/16/2015

Installation Operation Basic Maintenance



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 machine-readable form or any other form.

System IV Liquid to Liquid Heat Exchanger Instruction and Operation Manual

Table of Contents

PREFACE	
Compliance.....	3
WEEE/RoHS.....	3
After-sale Support.....	4
Warranty	4
Unpacking.....	4
Out of Box Failure.....	4
Feedback.....	4
SECTION I	
Safety	
Warnings.....	5
SECTION II	
General Information	
Description.....	6
Specifications	6
SECTION III	
Installation	
Site	7
Facility Water Requirements.....	7
Electrical Requirements.....	7
Voltage Selection	8
Plumbing Requirements	8
Fluids	9
Filling Requirements	10
Auto Refill Device (Optional)	10
Flow Control.....	10
Auto Restart.....	10
SECTION IV	
Operation	
Start Up.....	11
Temperature Controller	12
Tune Loop.....	13

SECTION V
Maintenance & Service

Service Contracts 14
Facility Water Strainer 14
Algae..... 14
Fault Interlock Contact..... 14
Configuration 15
Voltage Selection 15
Fuses 15
Phase Rotation 15
Pump Motor Overload Protector 16
Displaying Software Version 16

SECTION VI
Troubleshooting

Error Codes 17
Service Assistance 18
Technical Support 18

SECTION VII
Diagrams

Flow Diagram..... 19
Rear View 20

APPENDIX

Water Quality Standards and Recommendations..... 21

SECTION VIII
Warranty

..... 22

Preface

Compliance

Products tested and found to be in compliance with the requirements defined in the EMC standards defined by 89/336/EEC as well as Low Voltage Directive (LVD) 73/23/EEC can be identified by the CE Mark on the rear of the exchanger. The testing has demonstrated compliance with the following directives:

LVD, 73/23/EEC Complies with IEC/EN61010-1

EMC, 89/336/EEC IEC/EN61326-1

For any additional information, 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 both 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 or Service Contracts, contact our Service Department, see back cover. Before calling, *please* obtain the following information from the serial number label on the rear of the exchanger:

- BOM number _____
- Serial number _____

Warranty

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

Unpacking

Retain all cartons and packing material until the exchanger is 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 sellers 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 Thermo Scientific 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; seller to pay the cost of shipping. Customer must receive a Return Material Authorization (RMA) from Thermo Fisher Scientific prior to shipping the exchanger.

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 (see After-sale Support).

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.

Never operate the exchanger without cooling fluid in the reservoir.

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

Always empty the reservoir before moving the exchanger.

Never operate equipment with damaged line cords.

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.

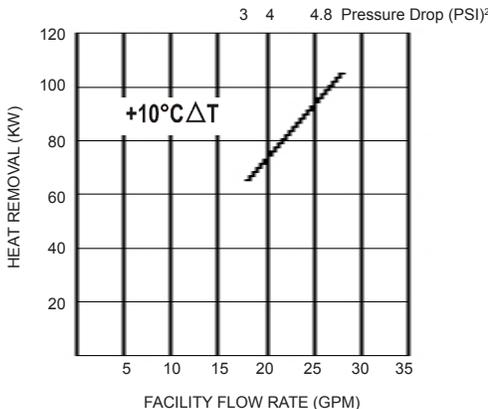
Section II General Information

Description

The Thermo Scientific System IV Liquid to Liquid Heat Exchanger is designed to remove heat from water-cooled instruments.

The System IV consists of a heat exchanger, recirculation pump, stainless steel reservoir and a microprocessor temperature controller.

Specifications

Temperature Range	+5°C to +40°C
Temperature Stability	±1.0°C
Heat Removal¹	
Cooling Capacity	100 kW / 341200 BTU
Pumping Capacity³	50 gpm @ 60 psi (189.3 lpm @ 4.1 bar) 41.5 gpm @ 42.1 psi (157 lpm @ 2.9 bar)
Reservoir Volume	
<i>Gallons</i>	6.5
<i>Liters</i>	24.6
Exchanger Dimensions⁴	
<i>(H x W x D)</i>	
<i>Inches</i>	33¾ x 23.0 x 27.0
<i>Centimeters</i>	85.7 x 58.4 x 68.6

Specifications obtained using water as the recirculating fluid and using water as a coolant on secondary loop, +25C 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 difference between the temperature of the cooling water supply (house water) and the cooling fluid as it leaves the System IV.
- Pressure drop indicates the minimum pressure differential pressure between the facility water inlet and outlet to achieve the corresponding facility water flow rate ($\text{Pressure}_{\text{inlet}} - \text{Pressure}_{\text{outlet}} = \text{Pressure}_{\text{drop}}$).
- If the pump trips or overloads it may be because the system pressure is too low or the pump flow is too high. Contact our Customer Support.
- Additional dimensions are shown on page 20.

Section III Installation

Site

Place the exchanger in a laboratory or clean industrial environment with easy access to a facility cooling water and a drain.

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 IV 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 meet these standards, the cooling capacity will be derated. The chart is based on a 10°C difference between the temperature of the cooling water supply and the process fluid entering the instrument being cooled.

As the heat load increases, the required flow rate of the cooling water supply increases. For example, if the heat load is 60 kilowatts, only 20 gallons per minute is required to remove the heat. However, if the heat load is 80 kilowatts, 26 gallons per minute 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 returning 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 to identify the specific electrical requirements.

Make sure the voltage of the power source meets the specified voltage, $\pm 10\%$.

Voltage	Frequency	Amperage	Line Cord Plug
230	60	14	L13-30
230	60	14	L15-30
380	50	6.9	None

Voltage Selection (60 Hertz Exchangers)

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

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

Plumbing Requirements

The plumbing connections are located on the rear of the exchanger and are labelled FACILITY WATER and RECIRCULATING WATER. FACILITY WATER connections are 1 inch nominal FPT. RECIRCULATING WATER connections are 1½ inch nominal FPT. See illustration on page 20.

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.

Connect the FACILITY WATER SUPPLY to the cooling water source.
Connect the FACILITY WATER RETURN to the drain.

Connect the RECIRCULATING WATER SUPPLY to the inlet of the instrument being cooled. Connect the RECIRCULATING WATER RETURN to the outlet of the instrument being cooled.

A basket strainer is supplied with the exchanger to protect it 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.

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 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.

If substantial lengths of cooling lines are required, they should be pre-filled with cooling fluid before connecting them to the exchanger.

Fluids



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

We recommend using distilled/deionized water with a 0.05 to 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

Remove the top access panel and locate the reservoir. Fill the reservoir with cooling fluid to within one inch of the top. Since the reservoir capacity is small compared to the volume of the circulating system, keep extra cooling fluid on hand until the circulating system is full.

Auto 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 solenoid 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.

NOTE: For the solenoid valve to close properly, the minimum supply flow rate should be 2.2 gallons per minute.

Connect the $\frac{3}{8}$ inch OD stainless steel barbed fitting on the solenoid valve to the make-up fluid source using $\frac{5}{16}$ or $\frac{3}{8}$ 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 “0” 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 exchanger will automatically restart when power is restored. This feature is enabled/disabled using the controller's Tune Loop, see page 13.

Section IV Operation

Start Up

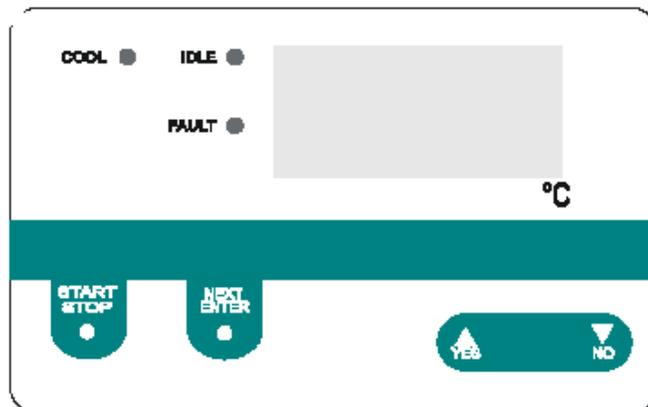
Before starting the exchanger, double check all electrical and plumbing connections and make sure the circulating system (the System IV, your application, and the tubing that connects them) has been properly filled with fluid.

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 the safe operating 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 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 RECIRCULATING 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 recirculation 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.

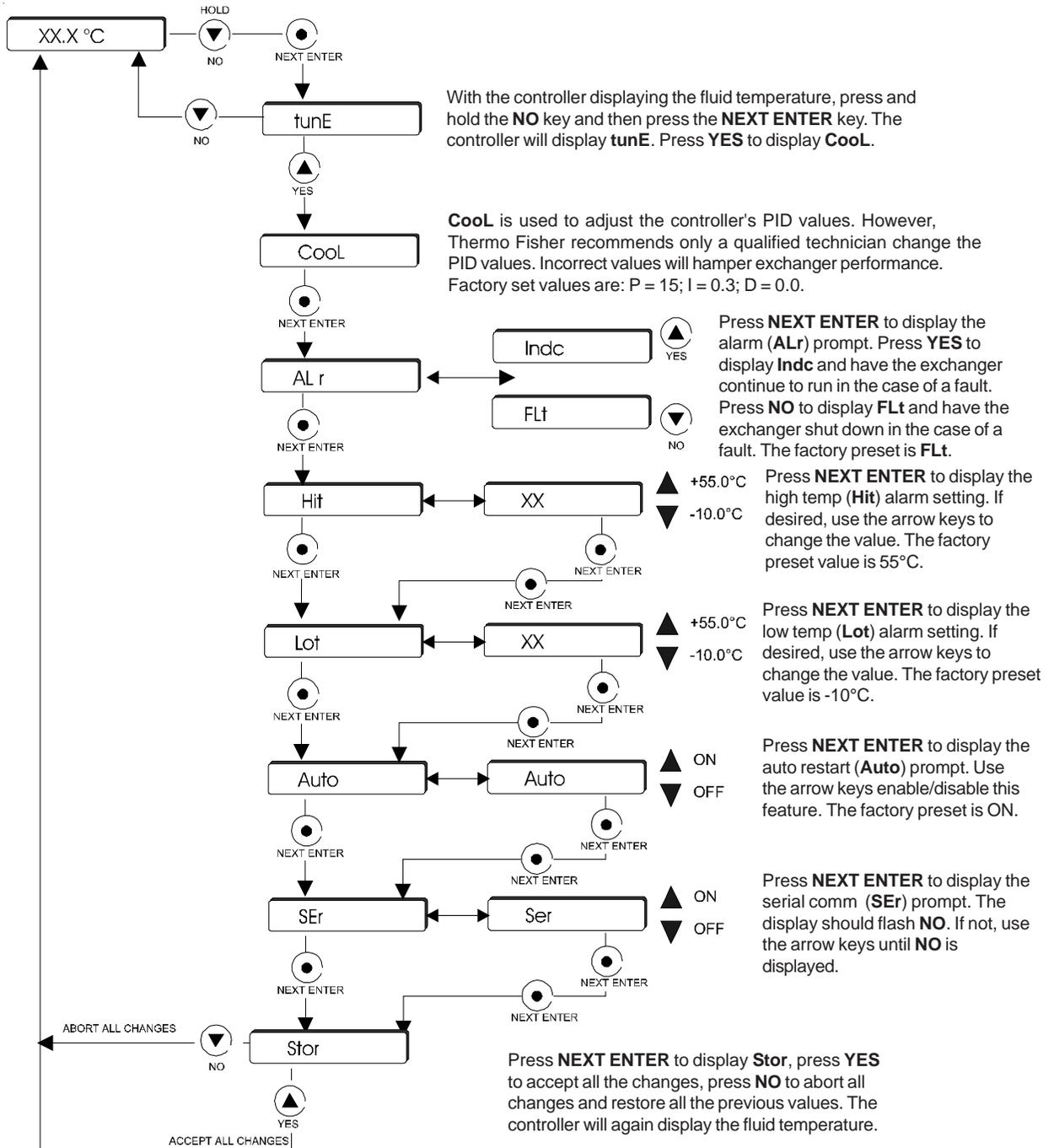
If the **NEXT ENTER** key is not depressed within one minute, the controller will time out and the new value will not be accepted. The controller will revert to the previous value.

NOTE: Error codes are addressed in Section VI, Troubleshooting.

TuneLoop

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.



Setup Loop

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).

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.

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.

Configuration

The exchanger top is secured to the cabinet by four ball stud retainers; one at each 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 may need to be removed. The access panels are secured using screws installed through the bottom of the cabinet.

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 as described in 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.

Fuses

The main fuses are 20 Amp RK5. The controller fuses are 1 Amp time delay. **Do not replace with an alternate amperage.**

Single Phase Exchangers

Remove the top cover from the cabinet.

The fuses are located on the right side of the control box.

Three Phase Exchangers

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 **OL** 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 microprocessor chip itself.

The following chart uses an example of software version 000550.93A

begin at the reservoir temperature display	20.3°C
Press and hold NO for at least 10 seconds.	0550
displays software version digits to left of decimal. Note the two leading zeros do not display.	
Press NEXT	93
displays software version digits to right of decimal.	
Press NEXT	1
displays software version revision letter (as its equivalent number - display cannot show letters. A=1, B=2, etc.)	
Press NEXT	0000
displays checksum - this can be disregarded	
Press NEXT	20.3°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 Er04 - Er15 persist, the problem is likely controller board failure. If Er16 is displayed the exchanger needs calibration, contact us for assistance.

Display	Indication
Er 04 - 13	Interrupt error
Er 14	Synchronous error
Er 15	Asynchronous error
Er 16	Bad calibration

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 Scientific customer service, see Preface.

Service Assistance

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

Part number

Serial number

Voltage of exchanger

Voltage of power source

Software Version

Technical Support

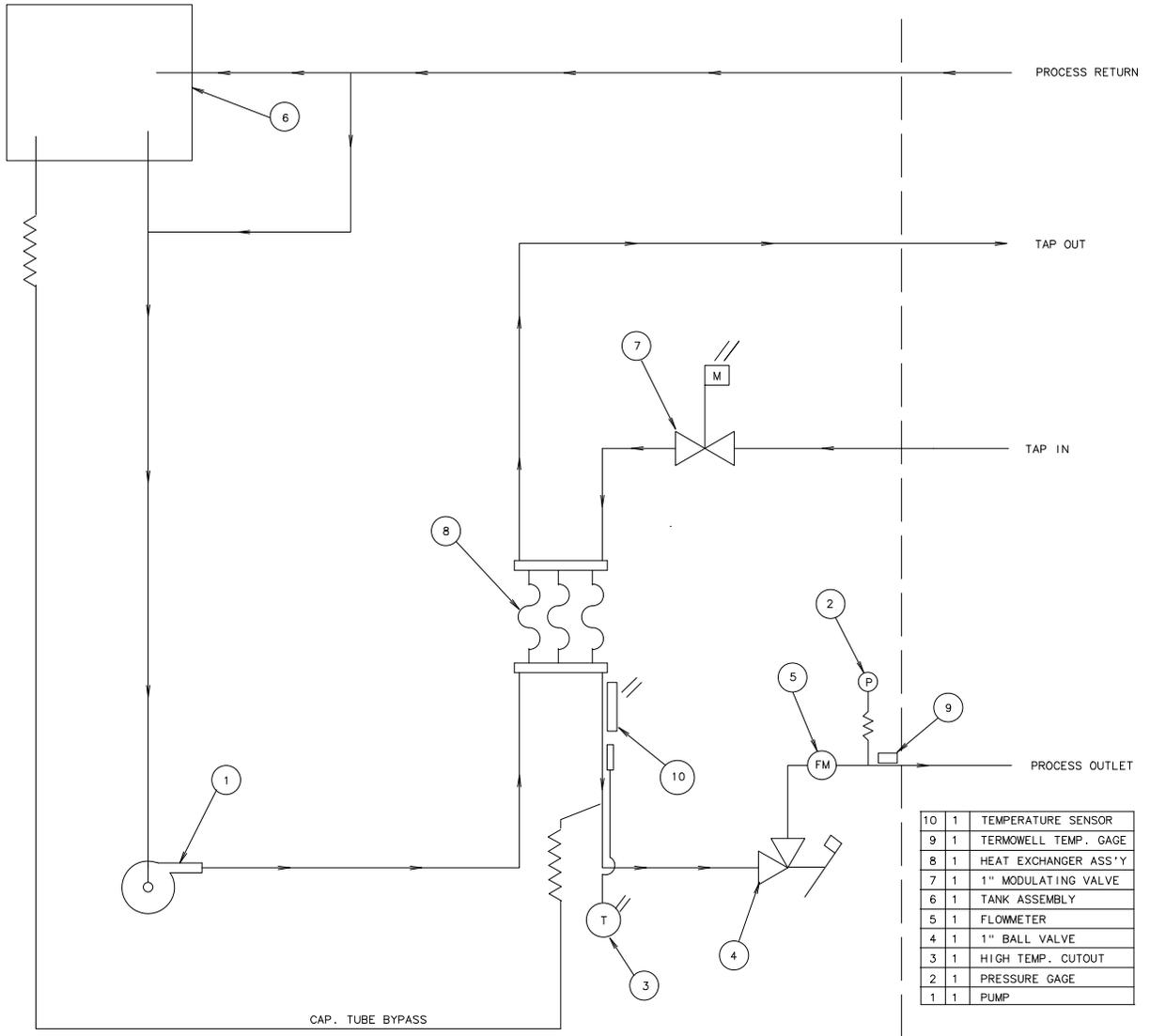
Our Service Department can provide you with a complete list of spare parts for your exchanger (see Preface, After-sale Support). Before calling, please obtain the following information:

Part number

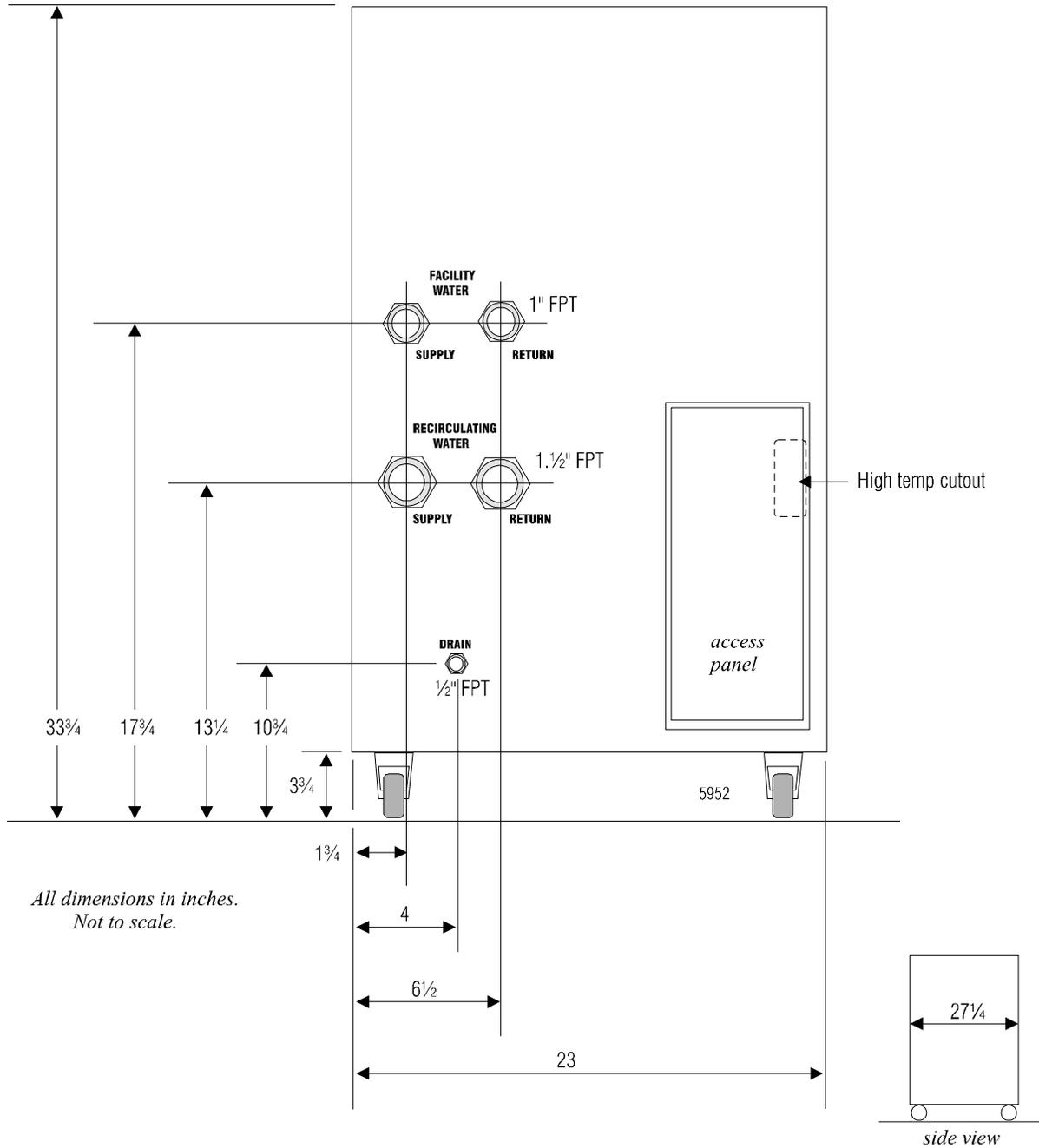
Serial number

Section VII Diagrams

Flow Diagram



Rear View



Appendix

Water Quality Standards and Recommendations

	Permissible (PPM)	Desirable (PPM)
Microbiologicals (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		
pH	6.5-8.5	7-8
Resistivity	0.01*	0.05-0.1*

* MOhmcm (Compensated to 25°C)

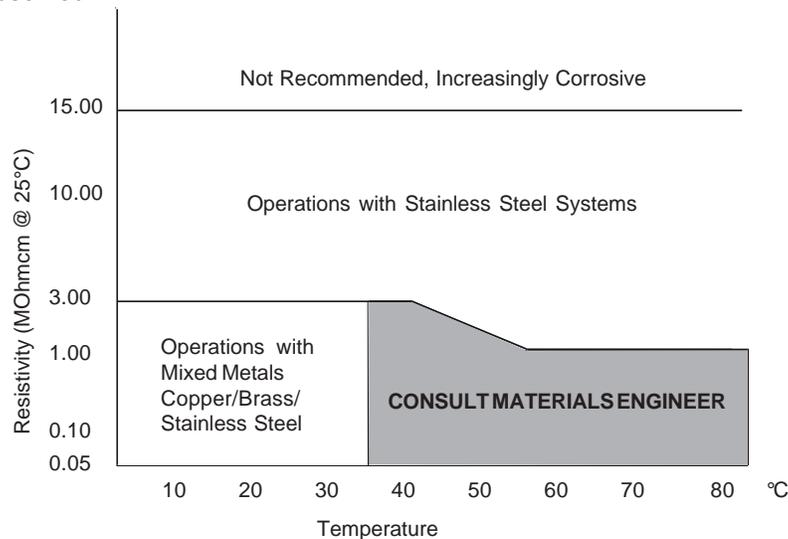
Unfavorably high total ionized solids (TIS) can accelerate the rate of galvanic corrosion. These contaminants can function as electrolytes which increase the potential for galvanic cell corrosion and lead to localized corrosion such as pitting. Eventually, the pitting could become so extensive that leaking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

High water hardness (Calcium and Manganese) 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 recommendations 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 rescission 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 Scientific ASSUMES 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