

CTL3000

Instruction Manual

Full Extractive Probe Controller Assembly

Part Number 111989-00

30May2008

© 2011 Thermo Fisher Scientific Inc. All rights reserved.

Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Thermo Fisher Scientific
Air Quality Instruments
27 Forge Parkway
Franklin, MA 02038
1-508-520-0430
www.thermoscientific.com/aqi

WEEE Compliance

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 Fisher Scientific products which may assist the detection of substances subject to the RoHS Directive are available at: www.thermoscientific.com/WEEERoHS.

Where to Get Help

Service is available from exclusive distributors worldwide. Contact one of the phone numbers below for product support and technical information or visit us on the web at www.thermoscientific.com/aqi.

1-866-282-0430 Toll Free

1-508-520-0430 International

We continue to support our customers with advanced online resources. Our Air Quality Instruments Online Library allows our customer's access to product documents and information on a constant basis.

Available 24-hours a day and seven-days a week, the online library provides quick access to information regardless of time zone or office hours.

To register for an account or log in, please visit www.thermoscientific.com/aqilibrary.

Claims for Damaged Shipments

A. The Thermo Fisher Scientific (TFS) CTL3000 probe controller assembly is shipped ready for operation. Immediate inspection of the TFS CTL3000 and probe controller assembly should follow upon receipt. Inventory of the container should be checked against the enclosed packing list. If there is a shortage of items, the operator should immediately contact TFS. If the contents are damaged, the carrier and TFS should be notified immediately.

B. The following documents are necessary to support claims:

1. Original freight bill and bill of lading.
2. Original invoice or photocopy of original invoice.
3. Copy of the packing list.
4. Photographs of damaged equipment and container.

NOTICE

Material contained in this manual is proprietary information of TFS. The TFS manual is to be used only for the installation, understanding, and servicing of this product. For further information or assistance contact:

TFS

All rights reserved. No part of this publication may be reproduced or stored in a retrieval system or transmitted in any form or by any means, electronic, mechanical photocopying, recording, or otherwise, without the prior written permission of TFS.

REVISION HISTORY			
Revision No.	Revision Made	Pages Affected	Revision Date
1.0	Initial release	All	May 30, 2008
2.0	Changed Company Name	All	July 10, 2009

1.0 PRODUCT DESCRIPTION 1-1

1.0	Product Description.....	1-1
1.1	Introduction	1-1
1.2	Probe Controller Assembly Description	1-1
1.3	Probe Controller Assembly Hardware	1-1
1.4	Specifications.....	1-2
2.0	THEORY OF OPERATION.....	2-1
2.1	General	2-1
2.1.1	Temperature Controllers	2-1
2.1.2	Solid State Relays.....	2-1
3.0	INSTALLATION AND OPERATION	3-1
3.1	Heated Filter and Heat-traced Temperature Controller Setup Procedure	3-1
3.2	General Installation.....	3-3
3.2.1	Power	3-3
3.2.2	Control and Data Lines	3-3
4.0	MAINTENANCE.....	4-1
4.1	General	4-1
4.2	Required Maintenance Equipment	4-1
5.0	ASSEMBLY REPAIR.....	5-1
5.1	Obtaining Replacement Parts	5-1
5.2	Spare Parts List	5-1
6.0	Appendix	6-1
6.1	XXXX7141, Rev 0, Sht 1 & 2, SOURCE PROBE CONTROLLER.....	6-1

LIST OF TABLES

3.1-1	Setup Menu Parameters.....	3-1
3.1-2	Operation Menu Parameters	3-2

1.0 PRODUCT DESCRIPTION

1.1 Introduction

The installation and operation manual provides instruction for basic installation, preventive maintenance, corrective maintenance, and trouble shooting procedures for the CTL3000 extractive probe controller assembly, which provides remote control of the PRO3000. This manual contains four sections:

- Section 1. Product Description: Hardware description, instrument operating parameters, and physical characteristics.
- Section 2. Theory of Operation: Complete functional description.
- Section 3. Installation and Operation: Instructions for installation and operation of the full extraction probe conditioning assembly.
- Section 4. Maintenance: Routine inspection, trouble shooting, corrective procedures, and repair/replacement for major assemblies.

1.2 Probe Controller Assembly Description

The TFS CTL3000 probe controller assembly is one part of a gas analyzing system. The system includes the PRO3000 probe assembly, the CTL3000 probe controller, one or more gas analyzers and a data acquisition system. The system operation and maintenance manual will describe the other parts of the system.

The CTL3000 probe controller assembly performs 3 functions:

- A) monitoring and control of the heated filter temperature in the probe conditioning assembly.
- B) monitoring and control of the heat-traced tubing umbilical temperature that transports the sample.
- C) connection of the extractive probe's alarm and control contact closures to the control system.

1.3 Probe Conditioning Assembly Hardware

The CTL3000 probe controller assembly enclosure mounts in a standard 19-inch rack and the total length of the enclosure is 15 inches. Additional inside connection clearance requires the rack to be at least 24 inches deep.

Connector #1 is used for connection to the extractive probe assembly. Connector #2 is for the connection to the rack controller. The back of the enclosure has a power entry module for a 120V AC cord connection.

The enclosure houses two temperature controllers and two solid state relays. These assemblies are discussed in Section 2.

1.4 Specifications

The TFS CTL3000probe controller assembly was designed to operate within the following specifications:

Power Requirements:	120V AC, 50/60 Hz, 30 Watts
Power Connection:	CSA/UL Approved power cord entry module
Operating Temperature Range:	-20°C (-4°F) to 50°C (122°F)
Weight:	18 lbs. (8.3 kg)

2.0 THEORY OF OPERATION

2.1 General

The CTL3000 probe controller assembly provides temperature control of the heated filter and tubing umbilical for the PRO400 extractive probe assembly and provides alarm contact closures to the customer's system.

2.1.1 Temperature Controllers

Two Watlow temperature controllers are located in the front of the CTL3000 probe controller assembly. Both of the controller readouts are in degrees Fahrenheit (°F) and can be changed to indicate degrees Celsius (°C) in the setup menu. The upper display is the actual temperature of the controlled device while the lower display is the temperature set point. On the face of each controller, located to the right of the temperature displays, are two LED's labeled "L1" and "L2." L1 indicates that power is being applied to a load. L2 indicates that the temperature has strayed past the alarm set point and an alarm will be sounding if a device is connected to the alarm output.

One of the controllers monitors and controls the probe heated filter temperature at a setpoint of 300°F by sensing the temperature of a resistance temperature device (RTD) located on the probe heated filter. The signal leads are connected directly to the temperature controller via connector CONN-1 on the rear of the chassis. The temperature controller outputs direct current to a solid state relay located in the controller. The solid state relay controls the heat cartridges mounted on the probe heated filter.

The same procedure is used for the heat-traced tubing umbilical with the exception of the solid state relay being located in the analyzer rack. The heat-traced tubing umbilical is temperature controlled to 300°F.

2.1.2 Solid State Relays

One solid state relay controls the band heater for the heated filter located in the probe conditioning assembly. A second solid state relay controls the purge valve located in the probe conditioning assembly.

3.0 INSTALLATION AND OPERATION

3.1 Heated Filter and Heat-traced Temperature Controller Setup Procedure


1. Remove the Watlow controllers from the front panel and check the dip switches on the rear of the controllers. Both switches should be to the OFF position.
2. Reinstall the temperature controllers and apply power to the probe controller assembly via the power switch located at the rear of the chassis.
3. Refer to Table 3.1-1 (below) for parameter set points for each controller. Enter the setup menu by pressing and holding the up and down arrows on the controller for three seconds. The controller will enter the setup menu when "LOC" is displayed on the lower window. To proceed to the next parameter, press the  (advance) key. *NOTE: all outputs will be switched off while in the setup menu.*

Table 3.1-1 Setup Menu Parameters

Parameter	Function	Heated Filter Settings	Heat Trace Settings
LOC	Lockout Security Level	0	0
In	Input Type	rtd	rtd
C_F	Centigrade/Fahrenheit	F	F
dEC	Decimal Point Placement	0	0
rL	Low Temperature Range	32	32
rH	High Temperature Range	572	400
Ot1	Output Type #1	Ht	Ht
HSC	Hysteresis Control	3.0	3.0
Ot2	Output Type #2	dEA	DEA
HSA	Alarm Hysteresis (Dead Band)	2	.2
LAt	Alarm Latching	NLA	NLA
rtd	RTD curve selection	din	din
rP	Ramping (not used)	OFF	OFF
PL	Power Limiting	100	100
dSP	Display Windows	NOr	Nor



4. Please refer to Table 3.1-2, below, for parameter set points for the controller. Enter the operation menu by pressing the  (advance) key. The controller is in the operation menu when “Pb1” is displayed on the lower window. To proceed to the next parameter, press the  (advance) key. These parameters will be automatically entered after the Auto-tune is complete..


Table 3.1-2 Operation Setup Menu

Parameter	Function	Heated Filter Settings	Heat Trace Settings
Pb1	Power Band #1	None*	None
rE1/dE1	Reset #1	None*	None*
Ra1/dE1	Rate #1	None*	None*
Ct1	Cycle Time #1	1.0	1.0
ALO	Low Alarm Setpoint	-10.0	-10.0
AHI	High Alarm Setpoint	+10.0	+10.0
CAL	Calibration Offset	0.0	0.0
Aut	AutoTune Function*	2 to Tune 0 for Off	2 to Tune 0 for Off

* Auto tune will be engaged when a value other than 0 is entered into the Aut parameter. When the Auto tune is complete, the value will automatically revert back to 0.

5. Verify that the sample pump is not energized. This will keep moisture from contaminating the system until temperature set points are reached.

7. Apply power to the extractive probe assembly by turning on the rotary power switch. Using a digital thermometer, check the temperature of the heated filter assembly. Ensure that the temperature is approaching the set point. When the temperature has stabilized, return to the probe controller assembly for the next step.

8. Ensure that both temperature controllers show a temperature within 10°F of the set point. Push the  (advance) key until “Aut” is on the display and use the “up” key to enter a “2.” The controller will begin an autotune and the heated filter temperature will cycle up and down until the controller has set autotune parameters. Perform the same function for the umbilical temperature controller.

9. Using a digital thermometer, verify the heated filter temperature is 300° F ± 5° F. Note that the thermal mass of the heated filter is large and may require 30-60 minutes to reach final temperature.

10. Verify that the sample pump is energized and running. Initiate a purge cycle on the extractive probe assembly.

3.2 General Installation

3.2.1 Power

A standard 120V AC, single-phase service is required to operate the CTL3000 probe controller assembly. Service is supplied using a #18 AWG power cord.

3.2.2 Control and Data Lines

Connect control lines from CONN-2 to the analyzer rack terminal strip using the specified shielded cable (refer to specific installation drawings for each application).

4.0 MAINTENANCE

4.1 General

The CTL3000 was designed to permit module replacement of failed assemblies. Other than replacement of failed assemblies or fuses, no maintenance is required. Failed controller boards may be returned to the factory for repair. To return a failed component to the factory, see Section 5.

4.2 Required Maintenance Equipment

To perform maintenance on the sample system, the following equipment is required.

- A. Normal hand tools.

5.0 RETURNING ASSEMBLIES FOR REPAIR

Should it become necessary to return any assembly, sub-assembly, or component for repair or replacement, contact the factory prior to shipment for specific information such as: return authorization number, shipping instructions, price, time to repair, etc. Also include pertinent facts describing the nature of the problem. Ship all components to the following address:

TFS

5.1 Obtaining Replacement Parts

The following information must be included in all purchase orders for parts:

- A. TFS Model and S/N of major assembly
- B. Part Number (found in parts tables)
- C. Description of part

5.2 Spare Parts List, CTL3000

TFS Part Number	Description	Recommended Quantity
22110019	CONTROLLER, TEMP., 93	1*
43005014	FUSE, 1/2 AMP DELAY MDL	1*
43003023	FUSE, 3 AMP DELAY MDL	1*
43002025	FUSE, 4 AMP DELAY MDL	1*

* THESE PARTS ARE RECOMMENDED AS A MINIMUM FROM THE TFS SERVICE DEPARTMENT TO AID IN ACQUIRING 98% UP TIME.

Printed in the U.S.A.

6.0 APPENDIX A

Drawing #	Rev	Sheet	Description
XXXX7141	0	1 of 2	SOURCE PROBE CONTROLLER
XXXX7141	0	2 of 2	SOURCE PROBE CONTROLLER

Warranty

Seller warrants that the Products will operate or perform substantially in conformance with Seller's published specifications and be free from defects in material and workmanship, when subjected to normal, proper and intended usage by properly trained personnel, for the period of time set forth in the product documentation, published specifications or package inserts. If a period of time is not specified in Seller's product documentation, published specifications or package inserts, the warranty period shall be one (1) year from the date of shipment to Buyer for equipment and ninety (90) days for all other products (the "Warranty Period"). Seller agrees during the Warranty Period, to repair or replace, at Seller's option, defective Products so as to cause the same to operate in substantial conformance with said published specifications; provided that (a) Buyer shall promptly notify Seller in writing upon the discovery of any defect, which notice shall include the product model and serial number (if applicable) and details of the warranty claim; (b) after Seller's review, Seller will provide Buyer with service data and/or a Return Material Authorization ("RMA"), which may include biohazard decontamination procedures and other product-specific handling instructions; and (c) then, if applicable, Buyer may return the defective Products to Seller with all costs prepaid by Buyer. Replacement parts may be new or refurbished, at the election of Seller. All replaced parts shall become the property of Seller. Shipment to Buyer of repaired or replacement Products shall be made in accordance with the Delivery provisions of the Seller's Terms and Conditions of Sale. Consumables, including but not limited to lamps, fuses, batteries, bulbs and other such expendable items, are expressly excluded from the warranty under this warranty.

Notwithstanding the foregoing, Products supplied by Seller that are obtained by Seller from an original manufacturer or third party supplier are not warranted by Seller, but Seller agrees to assign to Buyer any warranty rights in such Product that Seller may have from the original manufacturer or third party supplier, to the extent such assignment is allowed by such original manufacturer or third party supplier.

In no event shall Seller have any obligation to make repairs, replacements or corrections required, in whole or in part, as the result of (i) normal wear and tear, (ii) accident, disaster or event of force majeure, (iii) misuse, fault or negligence of or by Buyer, (iv) use of the Products in a manner for which

they were not designed, (v) causes external to the Products such as, but not limited to, power failure or electrical power surges, (vi) improper storage and handling of the Products or (vii) use of the Products in combination with equipment or software not supplied by Seller. If Seller determines that Products for which Buyer has requested warranty services are not covered by the warranty hereunder, Buyer shall pay or reimburse Seller for all costs of investigating and responding to such request at Seller's then prevailing time and materials rates. If Seller provides repair services or replacement parts that are not covered by the warranty provided in this warranty, Buyer shall pay Seller therefor at Seller's then prevailing time and materials rates. ANY INSTALLATION, MAINTENANCE, REPAIR, SERVICE, RELOCATION OR ALTERATION TO OR OF, OR OTHER TAMPERING WITH, THE PRODUCTS PERFORMED BY ANY PERSON OR ENTITY OTHER THAN SELLER WITHOUT SELLER'S PRIOR WRITTEN APPROVAL, OR ANY USE OF REPLACEMENT PARTS NOT SUPPLIED BY SELLER, SHALL IMMEDIATELY VOID AND CANCEL ALL WARRANTIES WITH RESPECT TO THE AFFECTED PRODUCTS.

THE OBLIGATIONS CREATED BY THIS WARRANTY STATEMENT TO REPAIR OR REPLACE A DEFECTIVE PRODUCT SHALL BE THE SOLE REMEDY OF BUYER IN THE EVENT OF A DEFECTIVE PRODUCT. EXCEPT AS EXPRESSLY PROVIDED IN THIS WARRANTY STATEMENT, SELLER DISCLAIMS ALL OTHER WARRANTIES, WHETHER EXPRESS OR IMPLIED, ORAL OR WRITTEN, WITH RESPECT TO THE PRODUCTS, INCLUDING WITHOUT LIMITATION ALL IMPLIED WARRANTIES OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE. SELLER DOES NOT WARRANT THAT THE PRODUCTS ARE ERROR-FREE OR WILL ACCOMPLISH ANY PARTICULAR RESULT.