



Ultra Low Temperature Freezers

Water Cooled Option

Installation and Operation

324003H02 • Revision A • September 2017

IMPORTANT Read this instruction manual. Failure to follow the instructions in this manual can result in damage to the unit, injury to operating personnel, and poor equipment performance.

CAUTION All internal adjustments and maintenance must be performed by qualified service personnel.

Material in this manual is for informational purposes only. The contents and the product it describes are subject to change without notice. Thermo Fisher Scientific makes no representations or warranties with respect to this manual. In no event shall Thermo be held liable for any damages, direct or incidental, arising from or related to the use of this manual.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

Contents

Introduction	1
Unpacking	2
Operating Requirements	3
Water Flow Rate.....	3
Water Pressure	3
Water Temperature Range	3
Water Quality	3
Water Drainage	3
Monitoring Equipment	3
Installation.....	4
Clearance.....	4
Water Connections	4
Operation.....	5
Maintenance	6
Cleaning the Water-Cooled Condenser	6
Replacing the Water Supply Line Strainer.....	6

Introduction

This manual describes installation and maintenance procedures for water cooling connections for Thermo Scientific Ultra-Low Temperature Freezers.



CAUTION: Be sure to read these instructions carefully before installing and using the water connections, and be sure to read the Ultra-Low Freezer Installation and Operation Manual before operating the freezer.

Use this product only in the way described in the product literature and in the user manuals. Before using it, verify that this product is suitable for the users intended application.

Unpacking

Upon unpacking, make sure that all specified water line fittings are included with the unit:

- 1/2" MNPT x 3/8" MNPT fitting (Quantity 2)
- 1/2" FNPT x 1/2" MBSPT adapter (Quantity 2)



Figure 1. 1/2" MNPT x 3/8" MNPT



Figure 2. 1/2" FNPT x 1/2" MBSPT

The 3/8" MNPT goes into the freezer at both the inlet and the outlet.

Operating Requirements

Water Flow Rate

The minimum water flow rate is 0.5 GPM (gallons per minute) if temperature of the water is less than 25°C and 1 GPM if temperature of the water is above 25°C.

Water Pressure

The water pressure should not exceed 90 psi or 620.5 kpa.

Note: This limit is not to be exceeded even though the components may be rated up to 150 psi.

Water Temperature Range

+12.0°C to +30.0°C (+53.6°F to +86°F)

Water Quality

Water must be free of particulates, which could cause a blockage or impair function of the regulating valve or heat exchanger. The pH level of water must be between 7.0 and 8.0. The optimum pH level is 7.4.

Use a suitable inline strainer in the inlet pipe to minimize particulates in the water supply. We recommend a strainer with a screen between 16 to 20 mesh, preferably stainless steel or brass. When the strainer is installed, water flow rate must still exceed the minimum water flow rates described in the **Water Flow Rate** section.

Water Drainage

If the water return line labeled PROCESS OUTLET is not connected to a closed loop chiller system, connect the return line water to a reservoir tank or drain. This will hold/dispose of the water from the unit to avoid flooding.

Monitoring Equipment

To verify operating requirements, you will need:

- A flowmeter capable of measuring up to at least 5.3 gallons (20 liters) per minute to measure flow rates at the water return line labeled PROCESS OUTLET.
- A pressure gauge ranging from 0 to 300 psig to measure pressure at the water inlet labeled PROCESS INLET.

Installation

Clearance

Before installing the water connections, be sure that the unit is positioned with an extra 2" (5 cm) of clearance in back to leave room for a hose.

This is in addition to the 6" (15 cm) back clearance recommended in the main user manual.

Water Connections

To install the water connections:

1. Uncap the PROCESS INLET/PROCESS OUTLET fittings on the freezer.



Figure 3. Water line connections on the freezer

2. Check fittings and remove any particles and/or debris.
3. Finger tighten the fittings, then turn 1/8 to 1/4 turn with an adjustable wrench.

Note: Thread sealant or Teflon tape is highly recommended for use on threaded fittings.

4. Check for leaks; if found, repeat the 1/8 turn to 1/4 turn until the leaks stop. **Do not overtighten.**



Figure 4. Fittings installed on freezer

5. Make sure the water line is clear from blockage and debris.
6. Install a strainer with a 16-20 mesh screen to the water-line PROCESS INLET.
7. Connect the incoming water supply line to the PROCESS INLET fitting.
8. Connect the water return/drain line to the PROCESS OUTLET fitting.
9. Adjust to specified operating standards (pressures, temperature).
10. Turn on the water flow to unit.
11. Check for any leakage.
12. Refer to the user manual for further startup instructions. **Make sure that you have read the main user manual carefully before starting up the freezer.**
13. Adjust the water flow rate. Note that water will not flow through the outlet until after the first stage compressor is running.

Note: Except for maintenance tasks described in the following section, do not turn on the freezer without water connected and flowing. You will not see water flowing through PROCESS OUTLET until after the first stage compressor is running.

Operation

Refer to the main ultra-low freezer user manual for all details about operation, parameter settings, general maintenance, user access, and troubleshooting.

Maintenance

The maintenance tasks described in this section should only be performed by qualified personnel.

Cleaning the Water-Cooled Condenser

Cleaning solutions can be used, depending on the type of deposits or build-up to be removed. Do not use liquids that are corrosive to stainless steel or the brazing material (copper or nickel).

The condenser should be cleaned at least once a year, more frequently when the environmental conditions are relatively high in particulates.



CAUTION: Do not use products containing hydrochloric acid or muriatic acid.



CAUTION: Cleaning solutions can be irritating to exposed skin. Take precautions as recommended by the manufacturer.

To clean the condenser:

1. Remove product from cabinet.
2. Disconnect the unit from the water supply.
3. Drain the unit by leaving it turned on until all of the water is out of the system, then turning the unit off **immediately**.
4. Hook up the water lines, turn on the unit, and rinse with fresh water.
5. Drain the unit again and turn it off.
6. Fill with fresh water and start the unit with water flowing.
7. Add cleaning agent to the water supply line labeled "PROCESS INLET". Choose the solution and concentration depending on the amount of deposits or build-up.
8. Circulate the cleaning solution.
9. Drain the cleaning solution.
10. Add and circulate a passivating liquid (phosphoric acid 5% or oxalic acid) for corrosion inhibition of plate surfaces. Drain this liquid.
11. Rinse with clear water and drain.
12. Reconnect the water supply and fill the unit.
13. Return to service.

Replacing the Water Supply Line Strainer

Recommended frequency of replacement depends on the operating conditions. If the water line being fed is clean, then the strainer should be replaced approximately every three months.

WEEE Compliance

WEEE Compliance. This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. It is marked with the following symbol. Thermo Fisher Scientific has contracted with one or more recycling/disposal companies in each EU Member State, and this product should be disposed of or recycled through them. Further information on our compliance with these Directives, the recyclers in your country, and information on Thermo Scientific products which may assist the detection of substances subject to the RoHS Directive are available at www.thermofisher.com under Services & Support.

Great Britain



WEEE Konformität. Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2012/19/EU erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliedstaaten und dieses Produkt muss durch diese Firmen wiederverwertet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Scientific, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identifizieren, die unter diese RoHS-Anweisung fallen, finden Sie unter www.thermofisher.com unter Services & Support.

Deutschland



Conformità WEEE. Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2012/19/EU in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Fisher Scientific ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Fisher Scientific con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo Scientific che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito www.thermofisher.com in Servizi e Supporto.

Italia



Conformité WEEE. Ce produit doit être conforme à la directive européenne (2012/19/EU) des Déchets d'Équipements Électriques et Électroniques (DEEE). Il est marqué par le symbole suivant. Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Fisher Scientific qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur www.thermofisher.com sous Services et Assistance.

France





IF YOU NEED ASSISTANCE:

Thermo Fisher Scientific products are backed by a global technical support team ready to support your applications. We offer cold storage accessories, including remote alarms, temperature recorders, and validation services. Visit www.thermofisher.com/cold or call:

Countries	Sales	Services
North America	+1 866 984 3766	+1 800 438 4851
India	1800 22 8374, +91 22 6716 2200	+91 22 6716 2200
China	+800 810 5118, +400 650 5118	+8621 68654588
Japan	+81 3 5826 1616	+81 3 3816 3355
Australia	+61 39757 4300	1 300 735 292
Austria	+43 1 801 40 0	+43 1 801 40 0
Belgium	+32 53 73 42 41	+32 2 482 30 30
France	+33 2 2803 2180	+33 2 2803 2180
Germany	0800 1 536 376, +49 6184 90 6000	0800 1 536 376
Italy	+32 02 95059 552	+39 02 95059 552, 432 254 375
Netherlands	+31 76 579 55 55	+31 76 571 4440
Nordic/Baltic/CIS	+358 9 329 10200	+358 9 329 100
Russia	+7 812 703 4215	+7 812 703 4215
Spain/Portugal	+34 93 223 09 18	+34 93 223 09 18
Switzerland	+41 44 454 12 22	+41 44 454 12 12
UK/Ireland	+44 870 609 9203	+44 870 609 9203
New Zealand	+64 9 980 6700	+64 9 980 6700
Other Asian Countries	+852 2885 4613	+852 2885 4613
Countries not listed	+49 6184 90 6000	+49 6184 90 6000

Thermo Fisher Scientific Inc.
275 Aiken Road
Asheville, NC 28804
United States

Find out more at thermofisher.com/cold

