thermoscientific



RVT450

Refrigerated Vapor Trap

Installation and Operation

80302105 • Revision A • May 2018



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

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

DISCLAIMER Thermo Fisher Scientific Inc. does not accept any responsibility for any damage caused to its products by unauthorized personnel.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

© 2018 Thermo Fisher Scientific Inc. All rights reserved.

Contents

Introduction	1
Safety Precautions	2
Operating Standards	3
Product Specifications	3
Environmental Conditions	3
Installation	4
Contents	4
Preparing The Refrigerated Vapor Trap	4
Operation	6
Emptying The Glass Flask	6
Accessories	7
Warranty	8

Introduction

Thermo Scientific Refrigerated Vapor Traps are reliable, easyto-use, benchtop cold traps for solvent vapor condensation and collection. Mechanically refrigerated vapor traps replace dry ice/methanol traps, and are the simplest, most practical, and safest way to protect high-vacuum, rotary vane oil pumps in a vacuum system.

Solvents are removed from samples during vacuum evaporation with a SpeedVac[™] Concentrator and are collected in the CFC-Free Refrigerated Vapor Trap. Solvent vapors pass into a Glass Condensation Flask and condense on its walls, preventing the vapors from reaching the oil pump and causing degradation.



CAUTION: To assure safe operation and best results, read this manual in its entirety before operating this instrument. Improper operation can damage the trap or your vacuum pump.

Safety Precautions

In this manual, the following symbols and conventions are used:



This symbol when used alone indicates important operating instructions which reduce the risk of injury or poor performance of the unit.



CAUTION: This symbol, in the context of a CAUTION, indicates a potentially hazardous situation which if not avoided could result in minor to moderate injury or damage to the equipment.



WARNING: This symbol, in the context of a WARNING, indicates potentially hazardous situations which, if not avoided, could result in serious injury or death.



This symbol indicates situations where dangerous voltages exist and potential for electrical shock is present.



The snowflake symbol indicates extreme low temperatures and high risk of frostbite. Do not touch bare metal or samples with unprotected body parts.



This symbol indicates a need to use gloves during the indicated procedures. If performing decontamination procedures, use chemically resistant gloves.



Before installing, using or maintaining this product, please be sure to read the manual and product warning labels carefully. Failure to follow these instructions may cause the product to malfunction, which could result in injury or damage. Below are important safety precautions that apply to this product:



WARNING: Disconnect the unit from all power sources before cleaning, troubleshooting, or performing other maintenance on the product or its controls.



WARNING: Do not use this device in radioactive, highly reactive or explosive atmosphere.

Do not use this device to process any explosive, radioactive, highly reactive or explosive atmosphere creating substances.

Operating Standards

Product Specifications

Temperature	-50°C (approx.)
Capacity	4 liters
Refrigerant	CFC free
Dimensions (W x D x H)	14 in x 24 in x 12 in
	36 cm x 61 cm x 30 cm
Weight	71 lbs
	32 kg
Operative Power*	115 VAC; 60 Hz; 8 amps
	230 VAC; 50 Hz; 6 amps
Fuse	8 A, 250 VAC, Time-lag
	6 A, 250 VAC, Time-lag

*Dependent upon ambient temperature, line voltage fluctuation, and load capacity.

Environmental Conditions

Indoor use only, in the absence of hoarfrost, dew, percolating water, rain and solar radiation.

Maximum altitude	2000 meters above mean sea level
Ambient temperature range	17°C to 32°C
Humidity	20% to 80% non-condensing
Pollution degree	2

Main supply voltage fluctuations not to exceed $\pm 10\%$ of the nominal voltage.

Transient overvoltages according to Installation Categories II.

Installation

Contents

- 1. RVT450 Refrigerated Vapor Trap
- 2. GCF400 Glass Condensation Flask
- 3. FC400 Flask Cover
- 4. M60-0044-11 Elbow Fitting
- 5. M60-0044-14 Straight Fitting

A second Glass Condensation Flask, GCF400, may be ordered separately. It is recommended that spare Glass Condensation Flasks are readily available, so one can be in use while others are being emptied, cleaned, and dried. CryoCool[™] Heat Transfer Fluid, SCC1 (1L) or SCC5 (5L), must be ordered separately.

Receiving. Immediately call manufacturer or your distributor if shipping carton shows any visible sign of damage.

Unpacking. Open the shipping cartons. Carefully remove the instrument and accessories. Lift and carry with two people, holding securely underneath with both hands. Use proper lifting technique (lift with the legs, not the back) to avoid personal injury. Compare the contents with the packing list. If there is a discrepancy, call Thermo Scientific technical service.

Inspection. Inspect the unit and accessories for damage that may have occurred during shipment. Should there be any damage, report it to the carrier and contact Thermo Scientific immediately. Make sure the carrier inspects the damage and leaves an inspection report. Register any claims for shipping damage against the carrier or his agent. Save the shipping carton in the event a return is necessary. Call Thermo Scientific technical service for further assistance.

Site preparation.

- The RVT450 requires a stable surface at least 26 inches (66cm) deep that is clean, dry, level, and within 4 feet (1.2 meters) of a compatible electrical outlet, 115 VAC & 60 Hz product should be plugged into a circuit rated for at least 8 amps, 230 VAC & 50 Hz product should be plugged into a circuit rated for at least 6 amps.
- The trap draws high current when first switched on, therefore, other high-powered equipment, or equipment that will be affected by a momentary drop in power should not be placed on the same circuit as the vapor trap.
- The product can also be placed on a sturdy cart, such as the Convenience Cart (CC120/DX), When used with a SpeedVac[™] Concentrator, we recommend placing the

concentrator on the top shelf of the cart, and placing the RVT450 directly underneath on the lower shelf. This placement will aid in trapping solvents efficiently.



CAUTION: Be sure to leave at least 4 inch clearance on all sides of the unit. Verify that the unit is on a leveled and stable platform. If necessary, move the unit to a more suitable location.



WARNING: Before connecting the unit to an electrical outlet, make certain that voltage, frequency, and amperage match the requirements indicated on the product label, name plate of the instrument. Use sockets with a protective earth conductor and correct mains cable.

Note: Do not use any detachable power cord that is not adequately rated for the unit.

Preparing The Refrigerated Vapor Trap

Some accessories are sold separately and must be purchased prior to set-up.

You will need:

- CryoCool™ Heat Transfer Fluid, SCC1 or SCC5 (sold separately)
- Spout Fitting, 1/2" x 3/4", M60-0044-24 (sold separately)
- 1. Switch the trap to **OFF**. Connect the power cord to the receptacle on the right side of the instrument. Plug the vapor trap into an appropriate wall outlet.
- Pour 750 ml of CryoCool[™] into the stainless-steel trap chamber up to the line scribed into the chamber wall. When replenishing CryoCool[™], add fluid until the level reaches the scribed line. CryoCool[™] is a non-toxic, nonflammable, long-lasting, odorless fluid that provides heat transfer between the flask and refrigerated chamber.
- Carefully place a clean Glass Condensation Flask into the chamber. Press down. Verify that the final CryoCool[™] level is 10–15 mm below the rubber seal. If the level is low, carefully pour more CryoCool[™] into the chamber while holding the flask in place. Immediately wipe clean any CryoCool[™] that spills onto the rubber seal.

4. To secure the flask in the chamber, insert the white Insulating Cover. Be sure the cut-outs for Flask Cover and finger notches face upward (see diagram).



5. Align the circular groove of the Flask Cover with the mouth of the Glass Condensation Flask. Press firmly into place to make a good seal.

Operation

After preparing the unit, switch it **ON** using the power switch located on the front panel (light will illuminate once turned on).



CAUTION: Immediately verify by touch that the trap is drawing air through the vent on the right side. If you cannot feel the air suction, switch the trap **OFF** at once. Operating the trap without a working fan, or with the air flow blocked, will damage the refrigeration system.

It may take up to 30 minutes for the trap fluid to reach its operating temperature. When this occurs, you may begin drying operations as outlined in the instruction manual of your drying apparatus.

The RVT450 is designed to operate continuously and may be left on for extended periods. Be sure to regularly check the Glass Condensation Flask and empty it as required. For maximum efficiency, replace with clean trap when not more than half-full.



CAUTION: Do not switch **OFF** and **ON** with in a short period of time as this could cause pressure to build up within the system and lock the system that could trigger the fuse. After switching **OFF**, allow sufficient time (minimum 15 minutes) for the pressure to stabilize before switching the system **ON** again.

Emptying The Glass Flask

During system operation, solvent vapors from the SpeedVac[™] will collect in the glass flask. This vessel must be regularly emptied to keep the entire system operating at peak efficiency. If the vessel is not maintained, it may become so full that sample drying rate is adversely affected. Another consequence of a full vessel is that solvent vapors are more likely to contaminate and possibly damage an oil-sealed vacuum pump that may be used with the SpeedVac[™] system.

Empty the glass vessel before it is half-full. Often it is convenient and good practice to change the vessel at the end of the workday, no matter the depth of fill. For aggressive solvents, you may wish to remove the vessel at the end of each run for maximum protection of the system components.



CAUTION:

 Use gloves while handling Glass condensation flask to avoid pain, local frostbites due to extreme low temperature.

- 2. Handle fully filled Glass condensation flask with care to avoid risk of injury.
- 3. Use face masks if required while cleaning the flask for protection against toxic chemicals and bio hazards.

To remove the GCF400 for cleaning, bleed the system back to atmospheric pressure. Remove rubber Flask Cover from flask, leaving tubing attached to cover. Withdraw the flask partially from the chamber and allow CryoCool™ to drain briefly. Fully remove flask and insulating foam ring. Avoid thermal shock by placing the flask on several thicknesses of absorbent paper toweling and allow to come to room temperature. Insert a spare GCF400, which is clean and dry, into the chamber. Cover with insulating foam ring, and seal with rubber Flask Cover. Make sure rubber cover is seated firmly for a good vacuum seal. This easy system maintenance can be done in a matter of minutes; the Refrigerated Vapor Trap need not be shut off during this process. When the used flask has defrosted, dispose of contents in an environmentally responsible manner per all applicable laws. Clean and dry flask for next use.

Note: If the Refrigerated Vapor Trap is not needed for several weeks, you may wish to shut it off between uses. Before switching on again, always remove the used glass vessel and replace with a clean, dry trap. Check the condition of the CryoCool[™]; if a layer of water (atmospheric condensation) is visible under the CryoCool[™], remove with a pipette. Failure to follow these precautions may cause the glass vessel to break when the trap returns to operating temperature



Note: Collected solvents may be purified for reuse or disposed of safely according to applicable regulations. If working with radioactive samples, test condensate for presence of radioactivity. If found to be radioactive, dispose of as radioactive liquid according to all applicable regulations.



WARNING: Disconnect the unit from all power sources before cleaning, troubleshooting, or performing other maintenance on the product or its controls.

Routine checks:

- 1. Check all the hoses to ensure that they are secure.
- 2. Ensure all glass condensation flask in emptied, cleaned and checked for crack before every run.
- 3. Clean the cover, cover seal and panels with a soft lint free cloth with one of the following
 - Mild detergent solution
 - Diluted Methanol (50%)
 - Diluted Ethanol (50%)

Accessories

- GCF400 Glass Condensation Flask, 4L, with Insulating Cover
- SCC1 CryoCool™ Heat Transfer Fluid, 1L
- SCC5 CryoCool™ Heat Transfer Fluid, 5L
- FC400 Flask Cover (black rubber)
- 145-6012-00 Insulating Cover (white)
- M60-0044-024 Spout Fitting, 1/2" x 3/4"

Warranty

All Thermo Fisher Scientific products mentioned in this manual (excluding glassware) are warranted against defects in workmanship for one year after the date of delivery to the original purchaser. This warranty is limited to defective materials and workmanship and does not cover incidental or consequential damages.

Thermo Fisher Scientific will repair free of charge any apparatus covered by this warranty. If a new component fails to work, Thermo Fisher Scientific will replace it, absorb all charges, and continue the one-year warranty period. Warranty work is subject to our inspection of the unit. No instruments, equipment, or accessories will be accepted without a Return Material Authorization (RMA) number issued by Thermo. Costs of shipping the unit are not covered under warranty. The warranty obliges you to follow the precautions in this manual.

When returning apparatus that may contain hazardous material, you must pack and label them following U.S. Department of Transportation (DOT) regulations applying to transportation of hazardous materials. Your shipping documents must also meet DOT regulations. **All returned units must be decontaminated (free of radioactivity, biological, or chemical contamination).**

Use of this equipment in manners other than those specified in this manual may jeopardize personal safety. Under no circumstances shall Thermo Fisher Scientific be liable for damages due to the improper handling, abuse, or unauthorized repair of its products. Thermo Fisher Scientific assumes no liability, express or implied, for use of this equipment.

Visit us online to register your warranty: www.thermoscientific.com/labwarranty

WEEE Compliance

WEEE Compliance. Products with this symbol are required to comply with the European Union's Great Britain Waste Electrical & Electronic Equipment (WEEE) Directive 2012/19/EU. 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/WEEERoHS under Services & Support.

WEEE Konformittät. Produkte mit diesem Symbol müssen der EU-Richtlinie für Elektro- und Deutschland Elektronik-Altgeräte (WEEE) 2012/19/EU entsprechen. Thermo Fisher Scientific hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen widerverwetet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Scientific, dieVerwerter und Hinweise die Ihnen nützlich sein können, die Thermo Fisher Scientific Produkte zu identizfizieren, die unter diese RoHS. Anweisungfallen, finden Sie unter www.thermofisher.com/WEEERoHS unter Services & Support.

Conformità WEEE. I prodotti con questo simbolo sono obbligatori per conformarsi alla Direttiva 2012/19/UE sui rifiuti di apparecchiature elettriche ed elettroniche (RAEE) dell'Unione Europea. Thermo Fischer 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/WEEERoHS in Servizi e Supporto.

Conformité WEEE. Les produits portant ce symbole doivent être conformes à la directive européenne relative aux déchets d'équipements électriques et électroniques (DEEE) 2012/19/UE. 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 le détection des substances sujettes à la directive RoHS sont disponibles sur www.thermofisher.com/WEEERoHS sous Services et Assistance.

Cumplimiento de la directiva RAEE. Los productos con este símbolo deben cumplir con la Directiva 2012/19/UE sobre equipos eléctricos y electrónicos usados (WEEE) de la Unión Europea. Thermo Fisher Scientific ha contratado a una o varias empresas de reciclado/disposición de residuos en cada estado miembro de la UE, y este producto debe reciclarse o desecharse a través de dichas empresas. Para obtener más información sobre nuestro cumplimiento con estas directivas, las empresas de reciclaje de su país, así como información sobre los productos Thermo Scientific que pueden ayudarle a detectar sustancias sujetas a la directiva RoHS, visite www.thermofisher.com/ WEEERoHS en la sección Servicios y Asistencia.







France



España





thermoscientific IF YOU NEED ASSISTANCE:

Thermo Fisher Scientific products are backed by a global technical support team ready to support your applications. Visit www.thermofisher.com/ or call:

Countries	Sales	Services
North America	+1 866 984 3766	(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/



©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. 80302105 0518