



Thermo Scientific Barnstead GenPure Ultrapure water system

Operating Instruction

50131265 Revision E June 2016

© 2016 Thermo Fisher Scientific Inc. All rights reserved.

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.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history:

For Research Use Only. Not for use in diagnostic procedures.

Preface

© 2016 Thermo Fisher Scientific Inc. All rights reserved.

These operating instructions are protected by copyright. Rights resulting thereof, particularly reprint, photomechanical or digital postprocessing or reproduction, even in part, are only allowed with the written consent of Thermo Electron LED GmbH.

This regulation does not apply to reproductions for in-plant use.

The contents of this operating instructions manual may change at any time and without any prior notice. Concerning translations into foreign languages, the English version of these operating instructions is binding.

Before you start to install and work with your ultrapure water system, please carefully read the information that is given in these operating instructions on how it is to be installed and operated.

This is particularly important as we, the manufacturer, cannot accept liability for any damage occurring as a result of incorrect operation of the system or from use of it for other than the specified purpose.

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.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Legal Information

NOTICE

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

Warranty

Thermo Electron LED GmbH warrants the operational safety and functions of the Thermo Scientific Barnstead Ultrapure Water Systems only under the condition that:

- the system is operated and serviced exclusively in accordance with its intended purpose and as described in these operating instructions,
- the system is not modified,
- only original spare parts, consumables and accessories that have been approved by Thermo Electron LED GmbH are used (third-party spares, consumables or accessories without Thermo Electron LED GmbH approval void the limited warranty),
- inspections and maintenance are performed at the specified intervals,
- an installation verification test is performed on commissioning the system for the first time and repeated after each inspection and repair activity.
The warranty is valid from the date of delivery of the system to the customer.
- The above mentioned warranty conditions are subject to the general terms and conditions of sale, in effect at the time of purchase, which apply as well.

Explanatory notes on the operating instructions



EU Mark of Conformity



CSA - admission



Indicates a situation which, if not avoided, could result in damage to equipment or property.



Indicates a potentially hazardous situation which, if not avoided, could result in death or serious injuries.



Indicates a hazardous situation which, if not avoided, will result in death or serious injuries.



General information! Particularly important notes are marked with this information sign.



Risk of electric shock! Electrical work on the system is only to be carried out by qualified personnel.



Protective conductor connection.
Connect the power supply to an electrical socket with a protective connection.



Indicates a situation where protected gloves or clothing is needed.



Indicates a situation in which protective goggles must be worn.



Indicates a situation in which breathing protection must be worn.

This information is valid for the system that is received.

For quick and correct service, please include the following information on all inquiries and replacement parts orders which relate to your system:

- The serial number (located on the right side of the system on the nameplate)
- The catalog number

Standards and Directives

The ultrapure water system complies with the following standards and directives:

- Low Voltage Directive 2014/35/EU
- EMC Directive 2014/30/EU
- ASTM D1193-6
- RoHs 2011/65/EU

Additionally, the ultrapure water system is in compliance with many other international standards, regulations and directives not listed here. Should you have any questions regarding compliance with national standards, regulations and directives applicable for your country, please contact your Thermo Fisher Scientific sales organization.

Contents

Preface	iii
Legal Informationiv
Warranty	iv
Explanatory notes on the operating instructions	v
Standards and Directivesvi
Chapter 1 Transport and packaging	5
Examination on receipt	6
Complaints	6
Packing for return shipment	6
Chapter 2 Safety precautions	7
Chapter 3 Extent of delivery	11
Extent of assembly kit	12
Available GenPure Systems	14
Chapter 4 Intended Use	15
Chapter 5 Technical specifications	17
Chapter 6 The installation area	21
Chapter 7 Installation	23
Start your system into operation	25
Wall mounting	27
Mounting the power pack (voltage supply)	29
Installation examples	31
Mounting example GenPure with Ion exchanger DI 1500 (option)	31
Chapter 8 Flow chart	33
Chapter 9 How the system functions	35
Chapter 10 Putting system into operation	37
Dispensing water from the dispensing valve	38
Venting the 0.2 µm final filter	38

Chapter 11	Operating elements	39
	Description of Display	40
	Flow chart of menu system control	41
Chapter 12	System control	43
	General information	44
	Operating modes	44
	Interval operating mode after switching on	44
	Non-stop mode	45
	Interval operation	45
	UV-Lamp	46
	OFF mode	46
	User menu	47
	Feedwater limiting value	47
	Ultrapure water limiting value	48
	UV-Lamp operating time and intensity	49
	Ultrapure cartridge operating hours counter	49
	Rinsing Procedure	50
	Disinfection Procedure	51
	Error history	52
	Print out of Data	52
	Entering a code number	53
	Code lock	53
	The OEM Menu	54
	Set the limiting value for temperature	54
	Set the rinsing time	55
	Change the disinfection time	55
	Set the interval pump time	55
	Set the interval rinse time	56
	Set the real-time clock	56
	Set the sending interval	56
	Language selection	57
	Switch units, conductivity/resistance	57
	Switch temperature compensation on/off	57
	Printer output	58
	Standard message	58
	Code message	58
	Error message	59
Chapter 13	Maintenance	61
	Maintenance intervals	62
	Change the ultrapure cartridge	63
	Disinfection	65
	Change the ultrafilter	68
	Structure of the UV-lamp	69
	Change the UV-lamp	71
	Change and autoclave the Final filter	75
	Autoclave the Final filter	76

Chapter 14	Waste disposal	77
Chapter 15	Trouble shooting	79
Chapter 16	Replacement parts	83
Chapter 17	Consumable materials	85
Chapter 18	Accessories	87
Chapter 19	Terminal assignments	89
Chapter 20	Maintenance records	91
Chapter 21	Contact Information Thermo Scientific	93
	Index	95

Transport and packaging

Contents

- “Examination on receipt” on page 6
- “Complaints” on page 6
- “Packing for return shipment” on page 6

 **DANGER**

Do not pull the plastic foil over your head. Risk of suffocation.
Use the plastic foil only for packaging.

Ultrapure water systems are carefully controlled and packed prior to dispatch, but damage could still possibly occur during transport. When the system is to be carried by hand, two people must always do this. Do not throw or tip the system.

Examination on receipt

- Check the completeness of the goods received against the delivery note.

NOTICE

Does the packaging show signs of damage? Inspect the system for damage.



Complaints

Should damage have occurred to the goods during transport:

- Immediately contact your delivery transport agency.
- Save the complete packaging, including the cardboard box, for a possible inspection of them and/or return shipment of the system.

Packing for return shipment

If possible, use the original box and packaging material.

When these are no longer available:

- Protect the system from shock by packing it in a suitable bag or sheet in a strong cardboard box.

NOTICE

The time limit for claims is 6 days from the time of receipt of the goods. The right to claim for damages ceases when this time has elapsed.

NOTICE

- Only a trained person is to be taken out the system out of operation.
- Prior to send back a operated system, empty the water and dry the system and take out the cartridges.
- Pack the filter cartridges into a bubble wrap and/ or packaging foam include it with the package of the ultrapure water system.

Safety precautions

NOTICE

Observe these safety precautions for your own safety.

CAUTION

Thermo Scientific Barnstead Ultrapure Water Systems are modern water purification systems intended solely for the treatment of potable water or water of ASTM Type II quality. The water it produces is not fit for drinking.

DANGER



Work may only be performed on the unit electronics when the unit has been switched off and when ESD protection is in place. Only specially trained personnel may work on the unit's electronics.

- Do not start to install and operate the system until you have read through the corresponding information given in these operating instructions.
- Lifting and carrying the ultrapure water system, e.g. to the installation location, should be carried out by two people. To lift it, each person takes hold of it under the base plate at two corners.
- The CE-mark is invalidated if constructional changes are made to the system, or if products of other manufacturers are installed in it.
- Protect the system from frost. The temperature in the area in which the system is installed is not to go below +2°C or above +40°C.
- Observe all appropriate rules and regulations, including the valid accident prevention regulations, which are applicable at the location where the system is installed.
- The feedwater pressure must be at least 0.1 bar and at max. 6 bar or 1.45 to 87 PSI. When the feedwater pressure is higher, install an external pressure reducer.
- A low pressure check valve is recommended to prevent back flow of feedwater from water system.
- A grounded 100-240V, 50/60Hz socket must be available (see “[Technical specifications](#)” on [page 17](#)).
- Access to the power supply cord and plug may never be restricted or obstructed.
- The installation area must have a drain at floor level with at least a nominal out diameter of 63mm or 12.48 inch (DN 50 pipe). Should no such drain be available it is recommend to install a water watcher (only for European specification). Otherwise the manufacturer will not accept liability for any possible water damage.

- If the system is to be at a rest for a longer time (e.g. long, holidays) switch the system off (unplug the mains plug) and shut off the feedwater line.
Allowing the system to run with the water feed line closed would result in damage to the pump. The manufacturer does not accept liability for such damage.
- Unplug the system from the power outlet for all Maintenance works on the system.
- The system must be subjected to rinsing and possibly also disinfection after longer rest periods. Follow the directions given in the section “Rinsing Procedure” on page 50.
- The surface or wall where the system is to be installed must have adequate load-bearing capacity (check the capacity and stability of the wall). The weight of the ultrapure water system is given under “Dimension and weight” on page 18.
- The surface on which the unit is installed must be level and stable - not to exceed a maximum of 2% deviation from evenness is recommended.
- When installing the ultrapure water system, always ensure that there is adequate space all around the unit (approx 30cm / 11.81 inch) to ensure that ease of use or replacement of materials (e.g. filter change, connection) is possible at all times.

WARNING

Never look directly into a switched-on UV-lamp, as UV-light endangers eyesight!

CAUTION

To avoid the risk of pinching, crushing cutting or electrical shock, never perform maintenance on the unit without its protective housing, or while it is in operation. Maintenance work on the system may only be performed by trained, authorized specialists.

- Visually inspect the system at regular intervals. Clean up any water or spills found around the system immediately



- Wear safety gloves when working with cleaning solution.
- If your skin should come into contact with a chlorine product, rinse it immediately with ample, fresh water.
- The unit or system components, may heat up as a result of a defect. It is recommended to always wear appropriate safety gloves to prevent skin damage or burns.
- Wear safety gloves when changing the UV-lamp, in order to prevent that your skin comes in contact with the UV-lamp glass.



- Wear safety goggles when working with cleaning solution.
- If your eyes come into contact with a chlorine product, rinse them immediately with ample, fresh water and immediately contact a physician.



- Check the UV-lamp before initial start.
 - If the UV-lamp is broken
 - wear directly a breathing protector (filter category FFP3) and replace the UV-Lamp
 - ventilate the room well.

NOTICE

The Hg content in the UV-lamp is so low so that no damage to the environment can arise.

- To avoid tripping ensure that the tubings do not lay over the floor.
- Apply the general rules of hygiene for laboratories when working with the system.
- Do not use oxidative cleaning agents when cleaning the system. They would cause damage to it.
- Proceed as follows when the system has a defect:
 - Switch the system off and unplug the system from outlet.
 - Shut off the feedwater supply.
 - Contact your local service organization.

Extent of delivery

Contents

- “Extent of assembly kit” on page 12
- “Available GenPure Systems” on page 14

Extent of assembly kit

Ultrapure cartridge
Catalog no.: 09.2005



Final filter 0.2 µm
Catalog no.: 09.1003

NOTICE

To increase the lifetime of the filter a sterilization at 120 °C for 30min is recommended. The procedure for the filter can be repeated up to 5 times.



Transformer-table power pack
Catalog no.:50149597



Universal Holder and Universal adapter
Catalog no.: 21.0007
Catalog no.: 21.0006



Connecting hose, 1.5m or 1.64 yards, straight/angle
Catalog no.: 18.0042



PE hose, 8x 1.2 m

Catalog no.: 18.0036



Connecting Cord (US) Catalog no.: 50132200

Connecting Cord (british) Catalog no.: 50132203

Connecting Cord (euro) Catalog no.: 50132215



Mounting parts for wall mounting the GenPure System:

- Plug 2 x S8

Catalog no.: 21.0035

- Screw hook 2 x 5.2 x 50mm or 5.2 x 1.97 inch

Catalog no.: 21.0057



Available GenPure Systems



50131211	GenPure standard	Basic system
50131235	GenPure UF	Basic system + ultrafiltration module
50131243	GenPure UV	Basic system + UV photooxidation
50131217	GenPure UV/UF	Basic system + UV photooxidation + ultrafiltration module
50131229	GenPure UV-TOC	Basic system + UV-photooxidation with TOC measurement
50131256	GenPure UV-TOC/UF	Basic system + UV-photooxidation with TOC measurement + ultrafiltration module

Intended Use

The Thermo Scientific Barnstead Ultrapure Water System is a laboratory system and is used for treatment of water. The system allows the purification of water into the water categories mentioned in the standards of ASTM 11.01 and ASTM 11.02.

The Thermo Scientific Barnstead Ultrapure Water System is designed to be installed and use in the following application areas:

- Laboratories for cell biological and biotechnological work with the safety levels L1, L2 and L3.
- Medical and microbiological laboratories according to DIN EN 12128.
- Laboratories in the central area of clinics and hospitals.

Unintended Use

The system must not be operated outside of the specifications as described in the operating manual. In particular, the system may not be used for production of drinking water and drugs manufacturing. The system must not be used as a medical device and outside of laboratories.

Technical specifications

NOTICE

Check at regular intervals the quality of your feedwater.

Demands the feedwater must fulfill

Source	Pretreated by reverse osmosis, ion exchange or distillation.
Clogging rate (SDI)	max. 1 for all versions. A 1 µm membrane prefilter is recommended for water not pretreated by reverse osmosis.
Feedwater resistance	> 0.5 MΩxcm
Free chlorine	max. 0.05 ppm
TOC	max. 50 ppb
Bacteria count	< 100 CFU/ml
Turbidity	< 1.0 NTU
Carbon dioxide (CO ₂)	max. 30 ppm
Silicate	max. 2 ppm
Particles	Filtration to 0.2 µm is recommended for protection of the internal filter / final filter.
Temperature	2 - 35°C
Pressure	0.1 - 6 bar or 1.45 to 87 PSI

Product water quality

		Standard	UV	UF	UV/UF	UV-TOC	UV-TOC/UF
Resistance (Reference temp. 25 °C)	MΩxcm	18.2	18.2	18.2	18.2	18.2	18.2
TOC	ppb	5 - 10	1 - 5	5 - 10	1 - 5	1 - 5	1 - 5
RNase	ng/ml	--	--	--	< 0.003	--	< 0.003
DNase	pg/ul				< 0.4		< 0.4
Bacteria	CFU/ml	< 0.1	< 0.1	< 0.1	< 0.1	< 0.1	< 0.1
Bacterial endotoxins	EU/ml	--	--	< 0.001*	< 0.001*	--	< 0.001*
Particles	µm/ml	< 0.2	< 0.2	< 0.2	< 0.2	< 0.2	< 0.2
Performance	l/min**	up to 2	up to 2	up to 2	up to 2	up to 2	up to 2

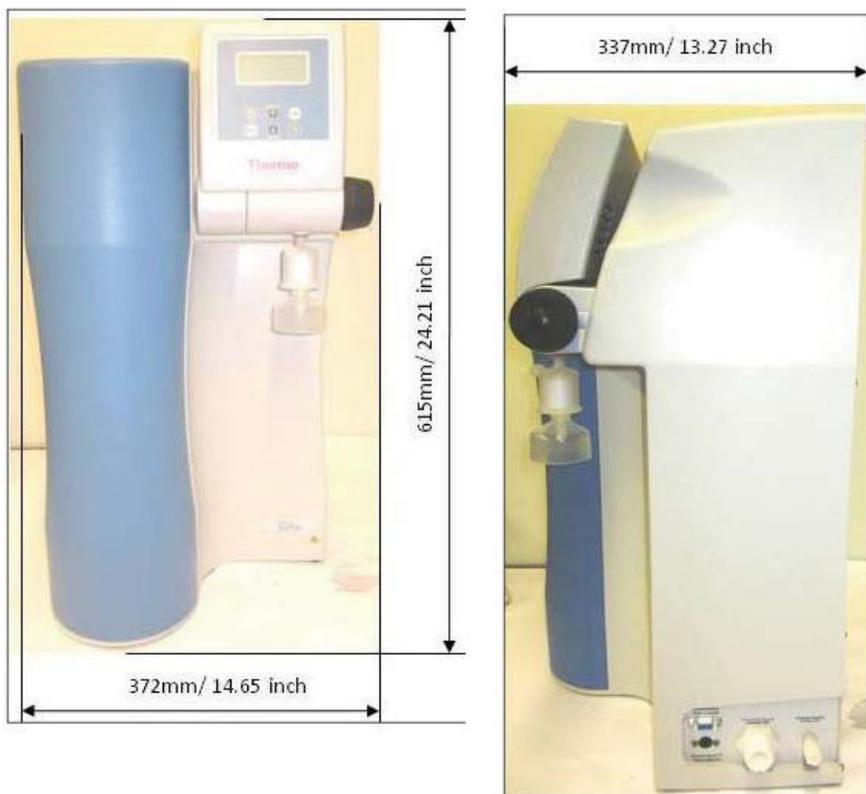
* Depends on the feedwater and disinfection

** Depends on the feedwater pressure

Dimension and weight		
Height	615 mm	24.21 inch
Width	372 mm	14.65 inch
Depth	337 mm	13.27 inch
Weight:		
GenPure Standard	22 kg	48.50 lbs (dry weight)
GenPure UF	23 kg	50.71 lbs (dry weight)
GenPure UV	24 kg	52.91 lbs (dry weight)
GenPure UV/UF	24 kg	52.91 lbs (dry weight)
GenPure UV-TOC	24 kg	52.91 lbs (dry weight)
GenPure UV-TOC/UF	25 kg	55.12 lbs (dry weight)

NOTICE

When the system is operating the system is by the amount of water about 3kg / 6.61 lbs heavier.



Cell constants of the measuring cells	
Feedwater conductivity	0.16 cm ⁻¹
Conductivity after UV oxidation	0.01 cm ⁻¹
Ultrapure water conductivity	0.01 cm ⁻¹
Connectors for water	
Feedwater	R 3/4"
Rinse water	8 mm OD hose
Ultrapure water	R 1/4"
Final filter outlet	8 - 10 mm OD hose
Electrical connections / external switched mode power supply	
Input voltage	AC 100 – 240VAC, 50 – 60 Hz, 2.0 A max
Output voltage	DC 24 V, 5.0 A max
System connection	DC 24 V, 80 W
Serial interface	RS 232
Protection Class	Class II (external SMPS certified as Class I)
Airborne sound emission	
Sound-pressure level	49 db(A)
Ambient conditions	
Operation area	Indoor rooms
Maximum altitude above sea level	Up to 2000 m
Temperature range during operation	min. +2°C, max +40°C, 80% rel. rH, non condensing
Temperature range storage	min. +2°C, max +60°C, 90% rel. rH, non condensing
Line-voltage variation	Not more than ± 10 % of the line voltage
Transient overvoltages	As usually occur in the supply network (overvoltage category II acc. to IEC 60364-4-443).
NOTICE	
The rated level of transient overvoltage is the withstand impulse voltage acc. to overvoltage category II of IEC 60364-4-443.	
Ventilation requirements	There are no special requirements with regard to ventilation.
Degree of pollution	2

Materials of parts which contact water	
Pressure reducer	NBR = Acrylnitril Butadien Rubber
Pump head	Nylon with glass fiber
UV-Lamp	High-purity synthetic quartz
UV Housing	Stainless steel
Ultrapure cartridge	PP = Polyethylene
UF Housing	PC = Polycarbonate
Rinsing solenoid valve	PA = Polyamide
Dispensing valve	PET = Polyethyleneterephthalate
Conductivity measuring cells	POM = Polyoxymethylen, stainless steel
Distributor block	POM = Polyoxymethylen
Connectors	POM = Polyoxymethylen
Hoses	PE = Polyethylene
O-Rings	EPDM = Ethylene Propylene Diene Rubber

The installation area

NOTICE

The operator is obliged to ensure, that the installation of the water purification unit and its operation are carried out in compliance with all national and international guidelines, applicable and valid for the place of installation. If necessary, measures to protect the drinking water have to be taken by installing appropriate components.

Take the following criteria into consideration when selecting the installation area:

Feedwater pressure, not below 0.1 bar (1.45 PSI) and not above 6 bar (87 PSI).

CAUTION

The feedwater pressure must never exceed 6 bar. If it is higher than this, install an additional external pressure reducer.

- Minimum air temperature + 2 °C.
- The surface on which the system is installed must be level and stable - not to exceed a maximum of 2% deviation from evenness is recommended.
- A smooth wall is required when the system is to be wall-mounted. Check the statics of the wall. It must have sufficient load-bearing capacity (for system weight, see “[Technical specifications](#)” on [page 17](#)).
- An atmospherically floor drain with a outer diameter of 63mm or 2.48 inch (DN 50 tube) shall be provided. When no floor drain is available, install a water watcher to protect against water damage (only available for EU). Otherwise the manufacturer will not accept liability for any possible water damage.
- Free run off to drain.

CAUTION

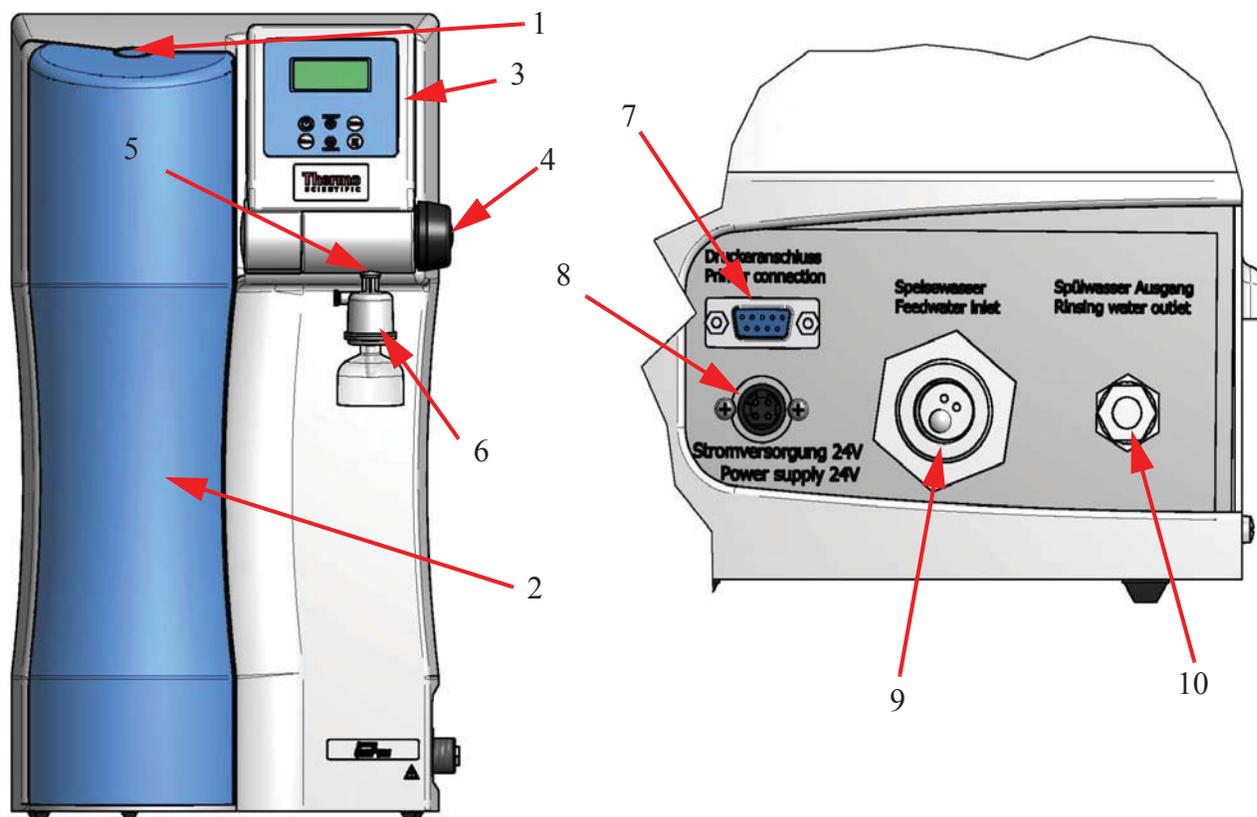
Unrestricted gravity flow to drain must be ensured!

- An electrical socket appropriate for the system (see “[Technical specifications](#)” on [page 17](#)).
- When installing the ultrapure water system, always ensure that there is adequate space all around the unit (approx 30cm / 11.81 inch) to ensure that ease of use or replacement of materials (e.g. filter change, connection) is possible at all times
- Easy access for operation and control of the system.
- Water pre treated such as DI, RO or distillation water connection with 3/4 NPT male thread and customer supplied shutoff valve.

Installation

Contents

- “Start your system into operation” on page 25
- “Wall mounting” on page 27
- “Mounting the power pack (voltage supply)” on page 29
- “Installation examples” on page 31

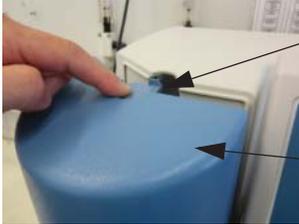
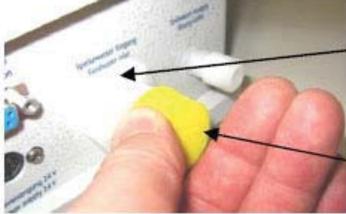


1. Push button for releasing the cartridge cover
2. Cartridge cover
3. Swivel Display
4. Rotary knob for ultrapure water
5. Connector for final filter R 1/4" female
6. Final filter 0.2µm
7. Optional printer connection
8. Power supply connector 24 V DC
9. Feedwater connector, R 3/4" male thread
10. Waste water connector 8mm o.d hose or 0.31" hose

Start your system into operation

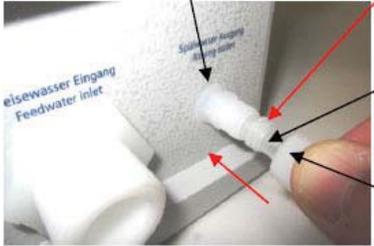
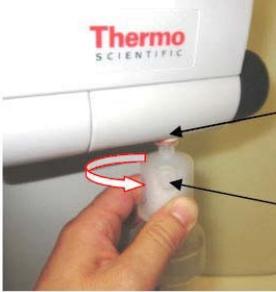
NOTICE

- Control after installation all tubings that the tubing have the right positions on the systems panel and there is no any leakage after open the feedwater supply tap.
- To avoid a book trip over the tubings, observe that the tubings are not lay over the floor.

Step	Action	Figure
1	Either place the system on the intended surface or hang it on a wall. For wall mounting using the included wall mounting hardware.	See under chapter “Wall mounting” on page 27”.
2	Release the cartridge cover by pressing the push button.	 <p>Push-button Cartridge cover</p>
3	Locate the ultrapure cartridge and fit the cartridge into the system.	
4	<p>Push each of the quick connectors onto the cartridge. You will know they are attached when an audible “click” is heard.</p> <p>Fit the cartridge cover on again.</p>	 <p>Quick connectors Ultrapure cartridge Inlet Outlet</p>
5	<p>Unscrew the yellow stopper from the feedwater connection of the system. Mount the connecting hose with 3/4” female thread nut which is supplied in the assembly kit to the feedwater connector of the system.</p> <p>The other end of the hose with 3/4” female thread you are mount it to an ion exchanger or an pre-treatment (see under chapter “Mounting example GenPure with Ion exchanger DI 1500 (option)” on page 31).</p>	 <p>Feedwater connecting Yellow stopper</p> 
	<p>CAUTION</p> <p>Only feedwater that has been pretreated by reverse osmosis, ion exchange or distillation is to be used!</p>	

7 Installation

Start your system into operation

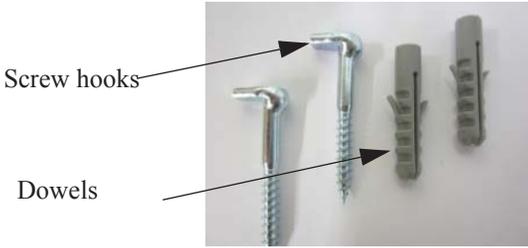
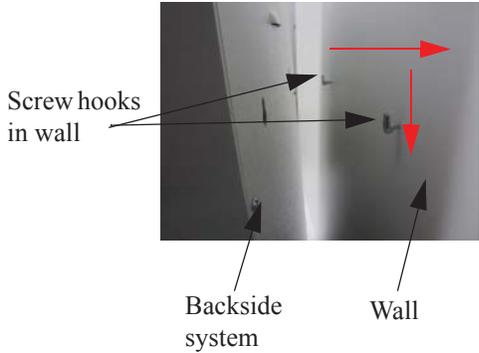
Step	Action	Figure
6	<p>Unscrew the fitting on the rinsing connector and pull the hose 8mm or 0.31 inch o.d through the fitting.</p> <p>Then push the white o-ring over the hose and screw the fitting back to the rinsing connector on the system.</p> <p>Make a gravity fall (pressure less) connection from the system (connector 10 to the floor drain). The drain to the sewer must be 1m or 1.07 yards above the rinsing water connector of the unit.</p>	<p>Rinsing water connector</p>  <p>White o-ring</p> <p>Hose 8 mm o.d</p> <p>Fitting</p>
7	<p>Screw the final filter supplied, counter-clockwise (see arrow in the picture) into the dispensing valve outlet (R 1/4" female thread).</p>	 <p>Connector for Autoclave final filter</p> <p>Final filter</p>
8	<p>Assemble the power pack and make the voltage connection to the GenPure system.</p>	<p>NOTICE</p> <p>See under chapter "Wall mounting" on page 27.</p>
9	<p>If applicable use the RS232 connector (7) to connect the optional data printer. Open the feedwater supply tap.</p> <p>CAUTION</p> <p>Only feedwater that has been pretreated by reverse osmosis, ion exchange or distillation is to be used!</p>	<p>Feedwater connecting kit</p> 

Wall mounting

NOTICE

You have the possibility to place your system onto a smooth surface or hang it on a wall. Before hanging the system onto a wall make sure that the wall can support the weight of the system once it's full of water.

Proceed as follows to hang your system onto a wall:

Step	Action	Figure
1	Draw with a pencil the distance from the holes to make the holes in the wall and then use a twist drill (8 mm or 5/16 inch) to make the two holes in the wall that are required as shown in the diagram.	See figure 1 holes for wall mounting
2	Plug the nylon S8 dowels (supplied in the assembly kit) in the holes. Screw the 5.2 x 50mm screw hooks into the dowels.	
3	<p>Lift the GenPure System and hang the back side of it onto the screw hooks.</p> <p>CAUTION</p> <p>Lifting and carrying the GenPure system should be completed by 2 people.</p>	

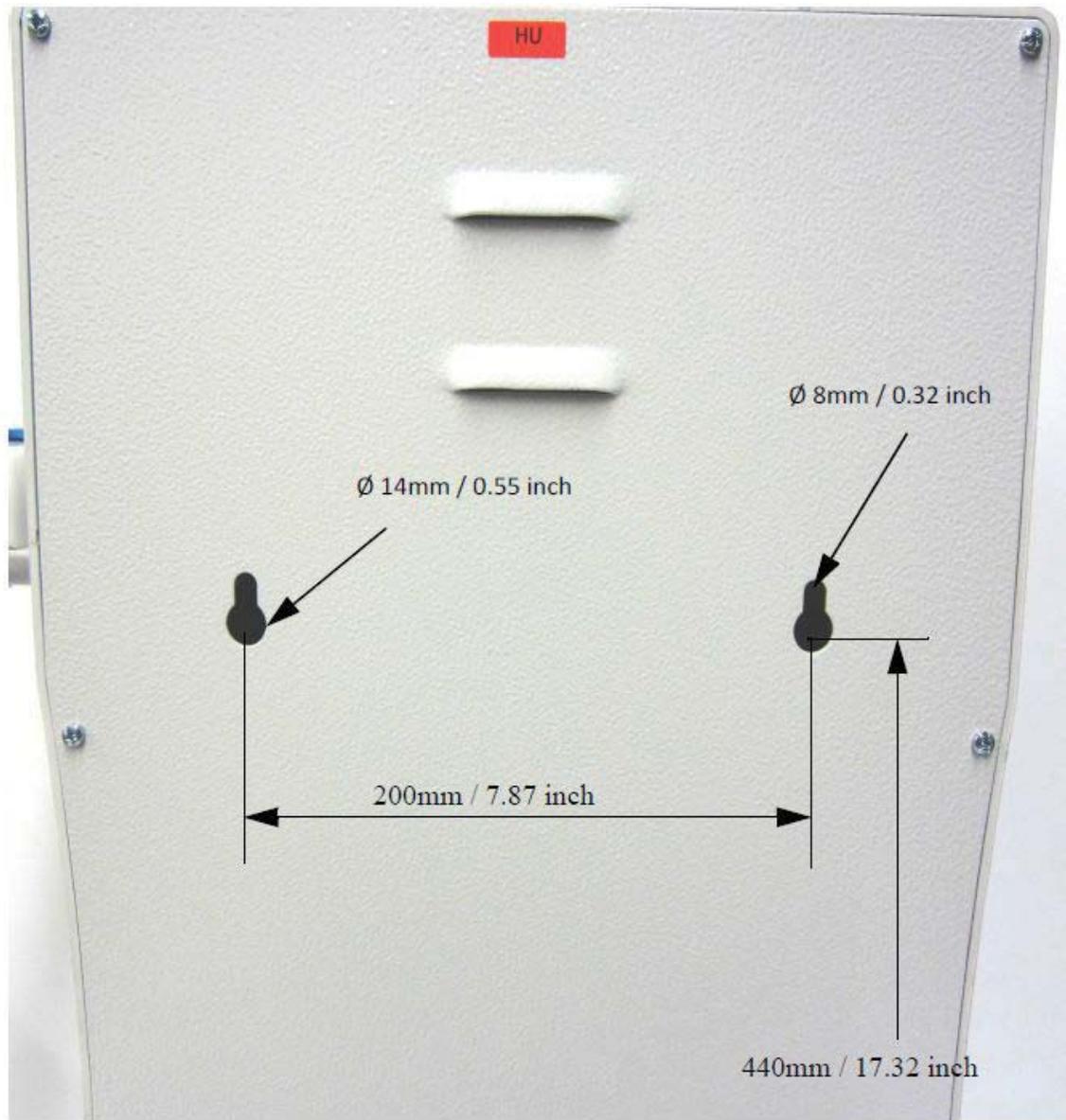


Figure 1. Holes for wall mounting

Mounting the power pack (voltage supply)

NOTICE

Whenever possible, mount the power pack on the wall to the left or right of the ultra pure water system where it is freely accessible and not come in contact with water for get wet.

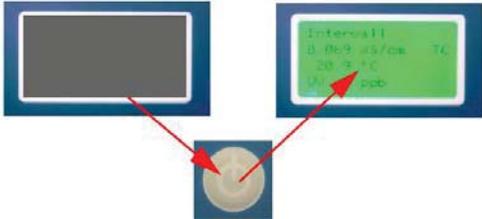
DANGER

Take caution to ensure that the suitable outlet and the power cable do not get wet and mount the power pack with dry hands.
Risk of an electrical shock.

Step	Action	Figure
1	<p>NOTICE</p> <p>Before beginning to work with the universal adapter and holder remove the protective foil from the backside of them.</p> <p>Stick the universal holder which is supplied in the assembly kit to the back of the power pack as shown in the figure next to this text.</p>	
2	<p>Stick the universal adapter to a smooth wall surface or screw it to the wall using the dowels and screws supplied in the assembly kit.</p>	
3	<p>When the universal holder and universal adapter have been fitted, hang the power pack in by pressing the power pack to the holder and then pull down (see red arrows).</p> <p>NOTICE</p> <p>The removable line cord must be shown to the bottom.</p>	
4	<p>Plug the connecting cable (appliance cable) in the power pack socket.</p> <p>DANGER</p> <p>Do not bring the power pack in contact with water. Risk of an electrical shock.</p>	

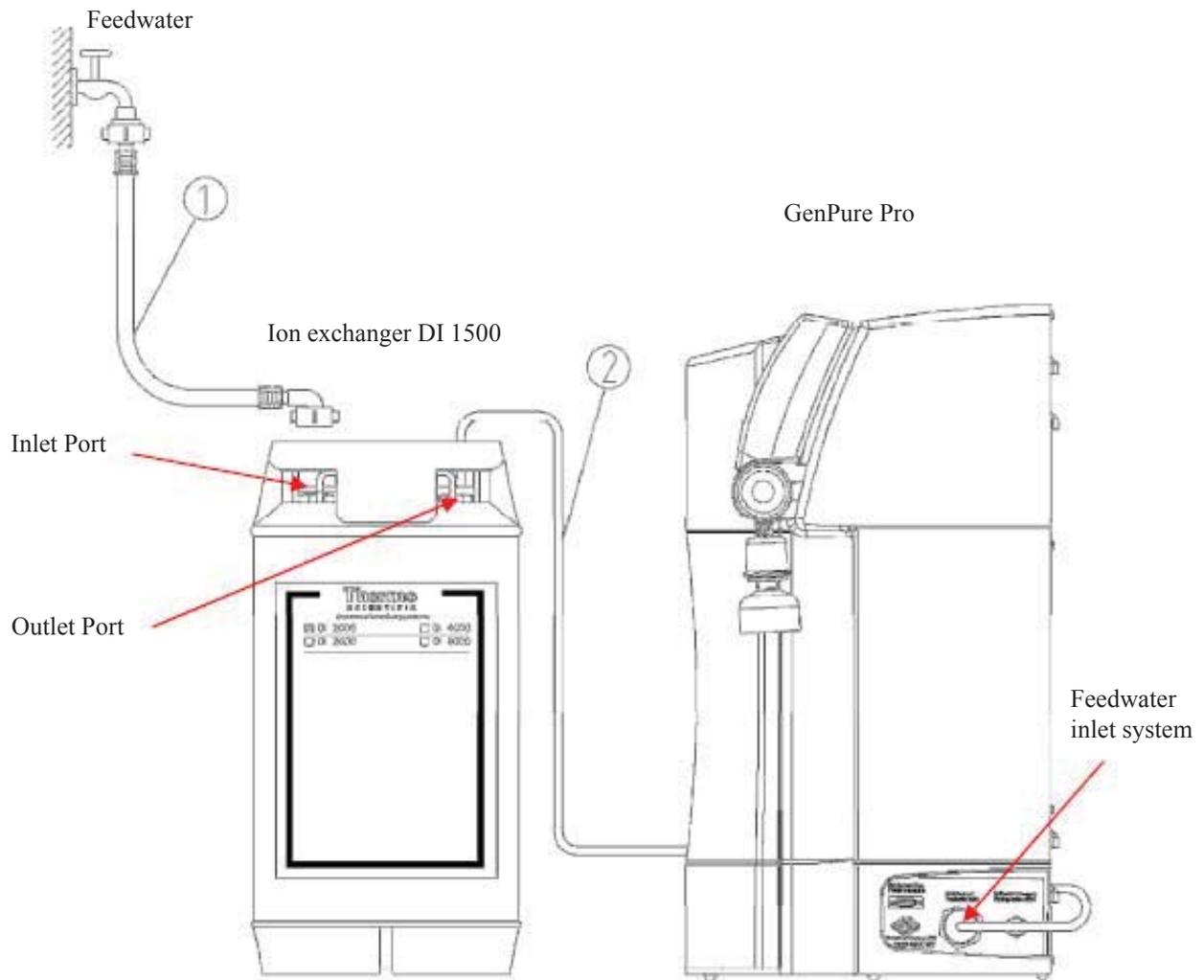
7 Installation

Mounting the power pack (voltage supply)

Step	Action	Figure
5	Connect the power pack to the ultrapure water system (24V 4-pin power supply connector, connector 3) and to an earthed 100 - 240V, 50/60Hz socket.	 <p>Power supply connector</p>
6	Switch the system on. The system is now ready for use.	

Installation examples

Mounting example GenPure with Ion exchanger DI 1500 (option)



Proceed as follows to connect an ion exchanger to the upstream side of the GenPure system:

Step	Action
1	Connect the hose which has a R3/4 female nut (1) from the raw water tap to the R3/4" input of the ion exchanger.
2	Make connection from the R3/4 output of the ion exchanger to the feedwater connector of the GenPure system by using the hose (2) that is contained in the assembly kit.

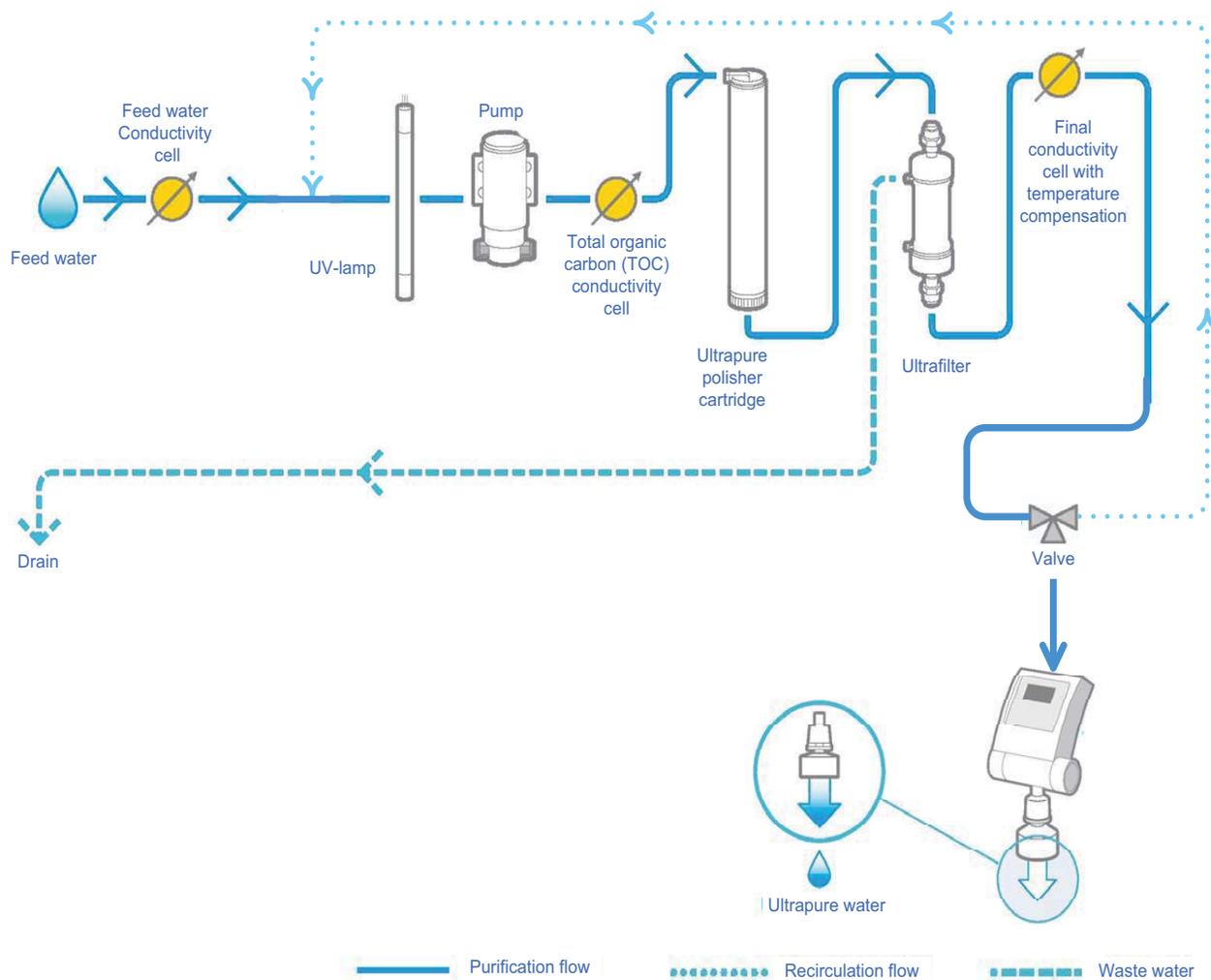
7 Installation

Installation examples

Flow chart

NOTICE

The following flow chart describes the GenPure system with full equipment (ultrafilter, UV-lamp, TOC measuring cell included). Depending on your GenPure system configuration the ultrafilter, UV-lamp or TOC measuring cell are inapplicable. The flow direction remains as described in the flow diagram.



How the system functions

NOTICE

System Function as applied in all GenPure systems

Tap water that has been pretreated upstream by reverse osmosis, ion exchange or distillation flows through a pressure reducer and into the ultrapure water system, where the conductivity is monitored. A pump directs this feedwater through UV-photooxidation (only possible in UV lamp equipped systems) and then through the ultrapure cartridge. From there the water flows through an ultrafiltration module (only possible in UF equipped systems). Then follows a permanent definition of conductivity measured by a special conductivity measuring cell equipped with temperature compensation. When ultrapure water is dispensed from the system, it flows through a final filter before reaching the point of use. During Interval operation, the water in the system is circulated in an internal circuit at regular intervals.

Systems with UV-TOC, UV-TOC/UF

Tap water that has been pretreated upstream by reverse osmosis, ion exchange or distillation passes through a pressure reducer and into the ultrapure water system, where the conductivity is monitored. A pump directs this feedwater through UV-photooxidation, which follows a conductivity measurement to determine the TOC value. Then follows an ultrapure cartridge and an ultrafiltration module (only with UV-TOC/UF), and the conductivity is then permanently measured by a special measuring cell (with temperature compensation). When ultrapure water is taken from the system, it flows through a final filter before reaching the dispensing outlet. During Interval operation, the water in the system is recirculated in an internal circuit at regular intervals.

The TOC value is calculated by taking the difference between the values measured by the measuring cells QIA300 and QI302. The measurement range is 0 - 30 ppb. When this range is exceeded, the number 99 is shown in the display instead of the measured value.

In Stand-by operation, “___” is shown.

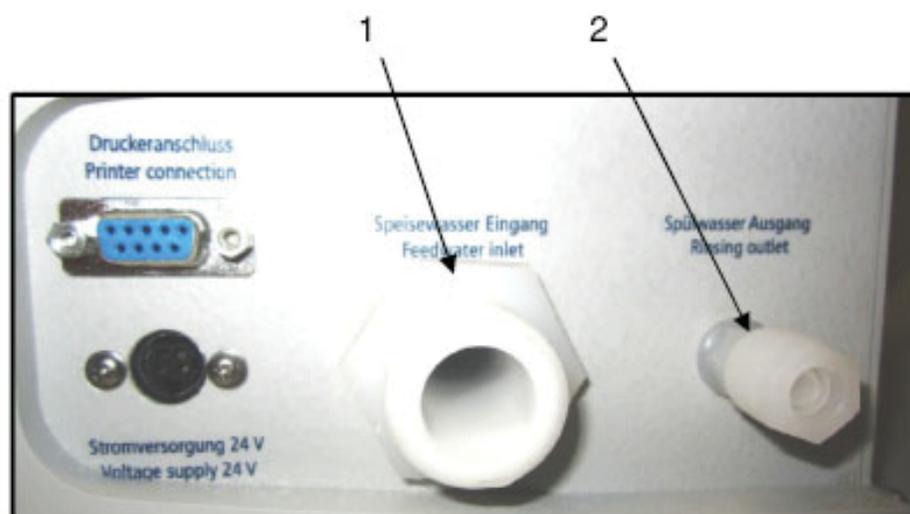
Putting system into operation

NOTICE

The system must have cooled down, or warmed up, to room temperature before being put into operation.

CAUTION

Check that all connections have been made as described above.



1. Feedwater connector R3/4" female thread
2. Rinsing water connector hose 8mm o.d or 0.31 inch.



Press this button to switch the system on. After a compulsory rinse, the system switches to the last used operating mode.

NOTICE

Vent the system by switching it to “Rinsing” three times in succession and, during this procedure, withdraw approximately 5 liters of water and discard it. The ultrapure water limiting value may be exceeded during this procedure.



Use the “NONSTOP” button to switch the system to the “Nonstop” operating mode“. This is the only mode which you can dispense water.



When the system successfully produces the ultrapure water quality that you require in “Nonstop” mode, press this button to return the system to the “Interval” mode.

Dispensing water from the dispensing valve

Step	Action	Figure
1	To dispense water from the dispensing valve push the dispense button on the handle once time. To stop dispense water from the dispensing valve push the dispensing button on the handle twice time.	

Venting the 0.2 μm final filter

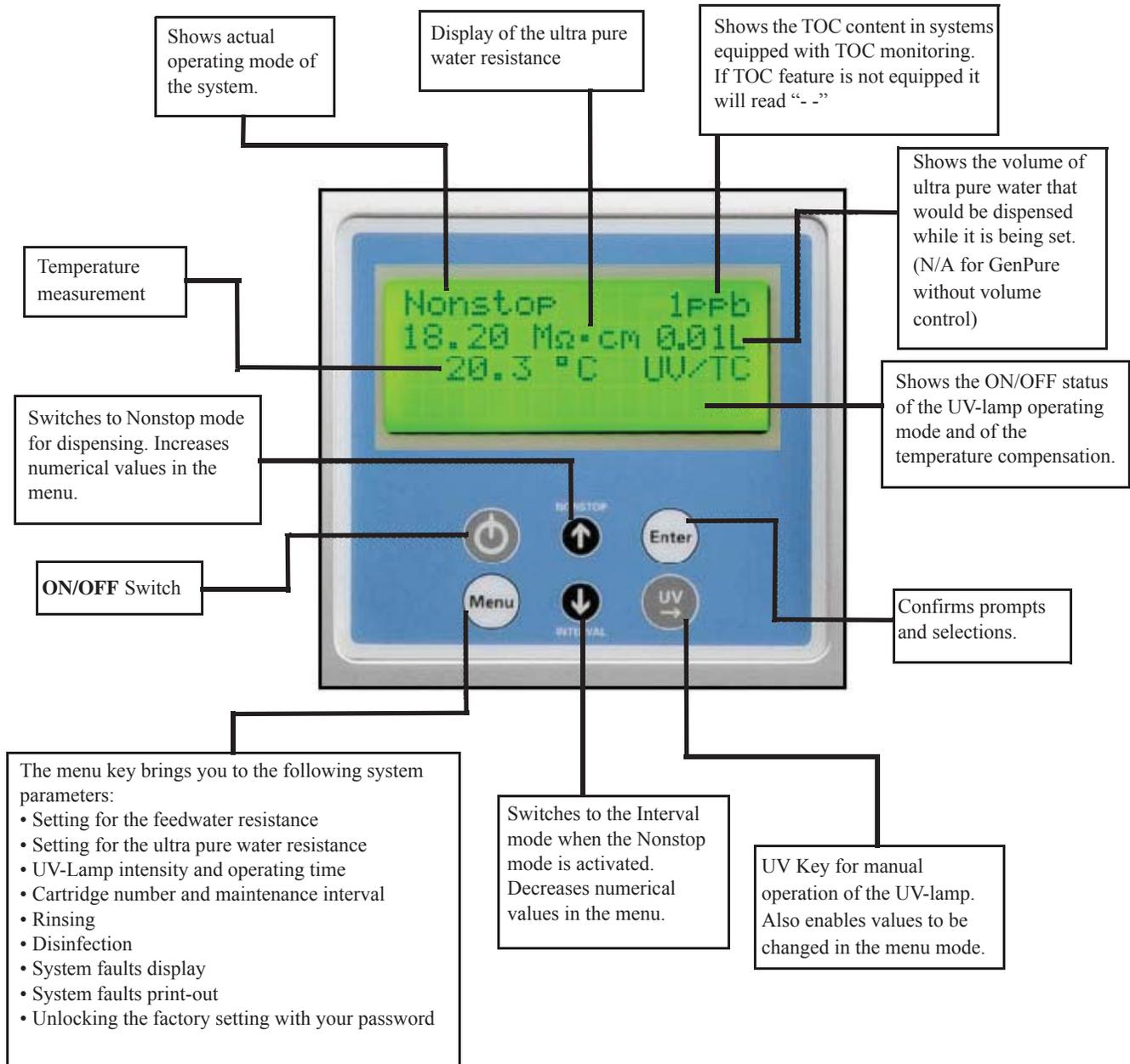
Step	Action	Figure
1	The first time you dispense pure water at the main dispenser through the 0.2 μm final filter, open the white knurled screw. NOTICE Do not close the knurled screw until pure water runs out of the opening at the knurled screw continuously. Rinse about 500 ml of water through the final filter.	

Operating elements

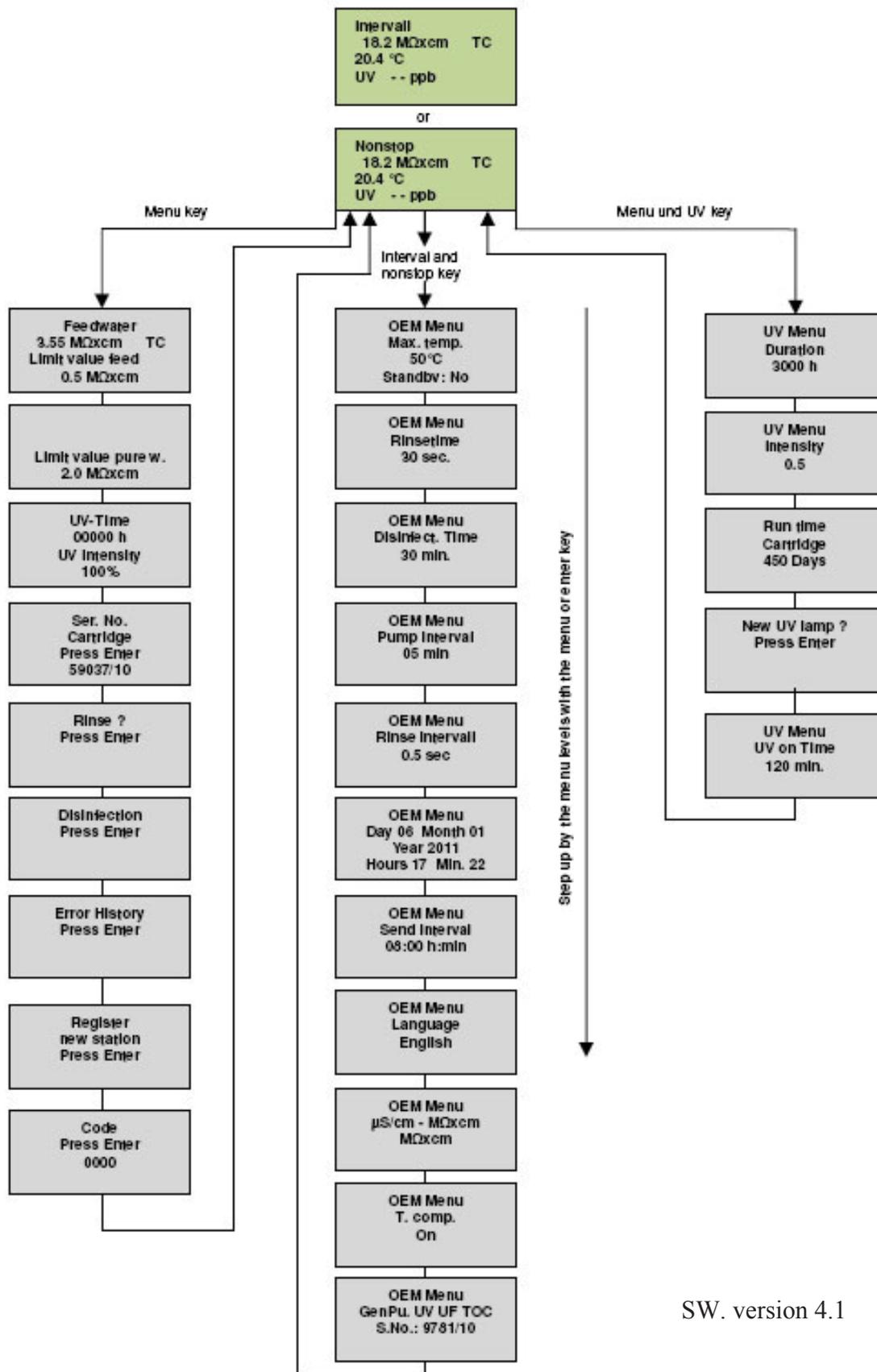
Contents

- “Description of Display” on page 40
- “Flow chart of menu system control” on page 41

Description of Display



Flow chart of menu system control



SW. version 4.1

System control

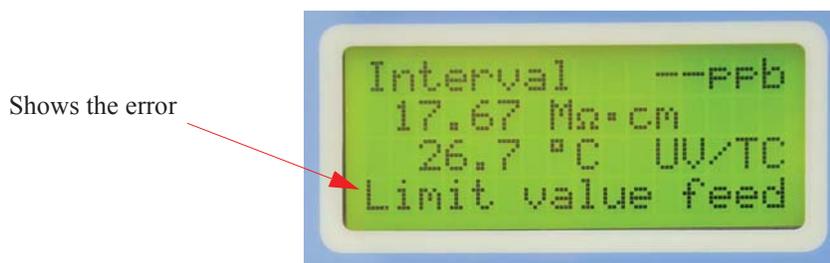
Contents

- “General information” on page 44
- “Operating modes” on page 44
- “User menu” on page 47
- “Code lock” on page 53
- “The OEM Menu” on page 54
- “Printer output” on page 58

General information

The software structure consists of five operating modes and four menus, which will be described in more detail in the following sections. Measured values are continually shown in the display and/or in the menus. The displayed TOC value is calculated from the difference in the ultrapure water measuring cell and TOC-measurement measuring cell values.

Should an error occur, the corresponding fault message is transmitted via the potential-free output and is shown in clear text in the 4th line of the display. In the case of several errors occurring at one time, they are alternately shown in the display.



Operating modes

Interval operating mode after switching on

Initially press the ON/OFF button. Then the display will show at first the system version, the system serial number and the software version number to display for 3 seconds. The system then automatically switches to the Interval operating mode (see “Interval operation” on page 45), whereby the green backlighting of the display is switched on and remains on until system control is switched off via the ON/OFF-button. The “UV” text message is displayed when the UV-lamp is switched on. The “TC” message is displayed when measured values are subject to temperature compensation. Further to these, the measured values for ultrapure water (measuring cell LF1) and temperature are also displayed. The displays of messages and measured values are independent of the operating mode.

The TOC value is not shown in Interval mode.
The display shows:



Non-stop mode

A press on the “nonstop” button switches the system to the non-stop mode. The non stop mode is the only mode in which water can be dispensed from the system. It is also the mode in which the system will continuously recirculate water with the system to keep the water ready for use. The circulation pump starts to run, the (UF) rinsing solenoid valve (V4) opens for the set “Intv.rinse time”. Non-stop operation is stopped automatically latest after 2 hours. Then the system operates in the “Interval”-Mode. The message *UV* is shown in the display when the UV-lamp is switched on. The UV lamp can only be switched on and off in this non-stop mode (see UV lamp). The TOC value is additionally shown in the display (TOC or *UV* only when applicable) whenever the UV-lamp is switched on for systems that have the TOC option.

The display shows:



Interval operation

The system is in the Interval mode when the system is switched on with the ON/OFF button. The Interval mode is used when the water system isn't needing to be in non-stop mode. This mode helps protect the system against bacteria growth as it will periodically recirculate water. Water can not be dispensed in this mode. The pump runs for the set interval pump time and the rinsing solenoid valve (V4) opens for the set “Interval rinse time”. When the interval pump time has expired, the pump is switched off until the end of the standstill time. The standstill time is given by the difference between half an hour and the interval pump time, so that the pump and the solenoid valve are actuated in a half-hourly rhythm. The TOC value is not shown in this operating mode.

The display shows:



UV-Lamp

A press on the UV-button results in showing the letters “UV”. However the UV-lamp is only switched on, however, when the system is in Nonstop operation. The UV-lamp is switched off at the end of Nonstop operation (settable). When Nonstop operation is manually ended by a press on the “Non stop” button, the UV-lamp is switched off after glowing for 0.5 hours. During the time that the UV-lamp is glowing. Furthermore the UV light intensity is monitored and is displayed in Menu (only applicable to systems with TOC monitoring). Should the limiting value for the UV-intensity (OEM menu / Menu) fall below a set value, the potential free output is set and the “UV Intensity” error message is displayed.

The operating time of the UV-lamp is recorded and the “*UV duration*” error message is brought to display when the limiting value set for this time is exceeded. TOC measurement is also carried out during the time that the UV-lamp is glowing only.

The display shows:



OFF mode

A second press on the ON/Off-button causes the display to go dark and all text output on the display to be extinguished. No outputs are now switched.

User menu

All measured values, operating times and limiting values which are relevant for the user can be set and read in this menu.

A press on the menu-button brings you to this menu. Each further press on the menu-button moves you further from one menu prompt to the next.

Settings can be changed with the arrow buttons. When you confirm a value by pressing on the Enter-button, you are taken to the next menu prompt. Settings can only be made when system control has been previously unlocked by entering a valid code number. (see “Code lock” on page 62)

To simplify changing settings, a press on the UV-button allows you to select a certain individual digit in the numerical value that you want to change. The arrow buttons can now be used to enter the new number from 0 to 9 at that position.

Feedwater limiting value

A single press on the menu-button allows the feedwater conductivity to be read or the limiting value of it to be changed. The fault message “Limit value feed” flashes in the 4th line of the display when the limiting conductivity value is exceeded.

Feedwater measuring range: 10 – 0.01 MΩ·cmΩ

Limiting value setting range: 0.1- 50.0 μS/cm

Basic setting: 0.2 μS/cm

Set the limiting value using the arrow buttons (see Settings with the arrow buttons).

With settings above 50 μS/cm, the limiting value is switched off and the word off appears in the display.

Press the Menu-button once then the display shows:



Ultrapure water limiting value

Two presses on the menu-key in this menu allow the fault display for the pure water limiting value and the pure water limiting value to be set. As soon as the fault display is switched on, the fault will be displayed both in Stand-by mode and in Production mode. When the fault display is switched off, the fault is only displayed in Production mode. The “Lim. val.pure w.” message is displayed when the limiting value is exceeded.

Ultrapure water measuring range:	0.1 MΩ·cm
Limiting value setting range:	0.055- 5.000 μS/cm
Basic setting:	0.2 μS/cm
Basic setting, fault suppression:	On

When a setting above 5.0 μS/cm is entered for the limiting value, the limiting value is switched off and the word “*Off*” appears in the display.

Press the Menu-button twice then the display shows:



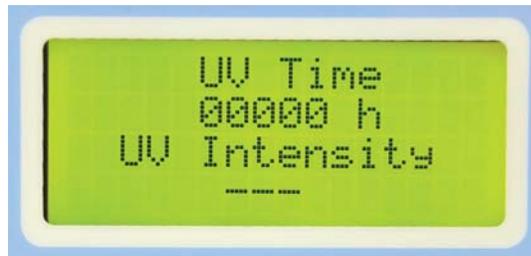
Ultrapure water measuring range:	10 – 0.01 MΩ·cmΩ
Limiting value setting range:	0.1- 50.0 μS/cm
Basic setting:	0.200 μS/cm

UV-Lamp operating time and intensity

In this menu the operation hours of the UV-lamp are indicated and the evaluation of UV-sensor input into the display under “*UV time*”.

The fault message “*UV duration*” is displayed when the maximum operating time has been reached. The UV-sensor measures the intensity of the UV-light, and this is displayed as a percentage value of the maximum value.

Press the menu-button 3 times then the display shows:



NOTICE

For more details see under section “Change the UV-lamp” on page 73.

Ultrapure cartridge operating hours counter

After fourth press on the menu-button the operating hours counter for the filter cartridge is set by input of a valid serial number.

Press the menu-button 4 times then the display shows:



NOTICE

For more details see under section “Change the ultrapure cartridge” on page 65.

Rinsing Procedure

A fifth press on the menu-button calls the question asking if rinsing is to be carried out. A press on the enter-button confirms this and triggers the rinsing procedure. The pump starts and the rinsing solenoid valve V4 opens for the rinsing time set in the OEM-menu.

The remaining rinsing time is shown in the display during rinsing.

Neither fault messages nor measured values are displayed during rinsing.

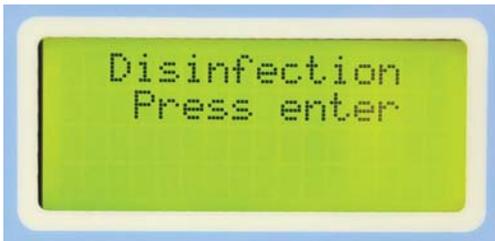
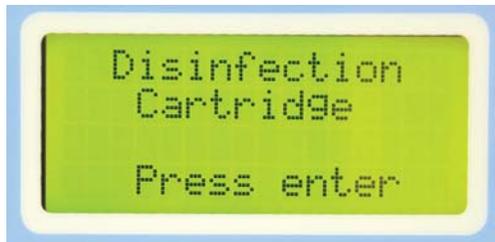
When the rinsing procedure is finished, the system returns to the last operating state (Interval or Nonstop).

Step	Action	Figure
1	Press the menu-button 5 times then the display shows:	
2	Confirm rinse by putting the enter button. The rinsing is started for 30 sec.	

Disinfection Procedure

A sixth press on the menu-button calls the question asking if a disinfection is to be carried out. A press on the enter-button confirms this, following which the demand “*Disinfection cartridge must be fitted*” is shown. When this has been fitted, a confirming press on the enter-button triggers the disinfection procedure. The pump starts for the full time set in the OEM-menu and, when the half of this time has elapsed, the rinsing solenoid valve opens and stays open until the disinfection procedure has finished. The demand “*New Filterset must be fitted*” is then displayed. When this has fitted, confirmation with the enter-button causes the system to return to the last operating state.

The remaining disinfection time is counted down and displayed during disinfection.

Step	Action	Figure
1	Press the menu-button 6 times then the display shows:	
2	Confirm disinfection by pushing the enter button. Change the filter cartridge with the disinfection cartridge (see under chapter “Disinfection” on page 67).	
3	Confirm with enter. The Disinfection is started for 30 min, indicating the remaining time.	

NOTICE

The completely process is described under section “Disinfection” on page 67.

Error history

Confirmation of this prompt with Enter allows the error storage to be looked through. Two errors, each with date and time, are shown in the display. Pressing the arrow buttons takes you successively through preceding or following errors. Press the menu-button to end the error display. This takes you to the next menu prompt.

Step	Action	Figure
1	Press the menu-button 7 times then the display shows:	
2	Confirm error history by putting the enter button. Now you can see two last saved errors with date and time. The error code can be requested at the local service organization.	

Print out of Data

In this menu, the current system data can be printed via a connected printer.

Press the menu-button 8 times then the display shows:



Entering a code number

To prevent unauthorized access to system control, settings can only be changed when a valid code number is entered and confirmed with Enter in this menu. Each code access is issued to the printer (RS 232) with date, time and code number. Valid codes are found in this manual in section “Code lock” on page 62.

Press the menu-button 9 times then the display shows:



NOTICE

You can assign the permissible code numbers listed in the Table on the following page to appropriate members of the staff etc.

When names have been entered, tear the page out and file it where it is safe from unauthorized viewing.

NOTICE

Press the menu-button 10 times and you are leave the User menu and the system is back to the last System system operation that you have choose.

Code lock

To prevent unauthorized access to system control settings, changes to these settings can only be carried out when a correct code number has been entered and confirmed with Enter.

In deviation to existing programs, control release can be given at three levels. Only the menu is released for changes at the first level. Both the menu and the OEM menu are released at the second level. All menus are released at the third level.

Code numbers:

No.	Menu	No.	Menu + OEM menu	No.	All levels
1	0150	4	0450	7	0750
2	0250	5	0550	8	0850
3	0350	6	0650	9	0950

Each access via the code is printed out by the printer (RS 232) complete with date, time and the code number used.

The OEM Menu

Basic settings and limiting values can be changed in this menu. To make such changes in the OEM-menu, the system must be unlocked first (see “Code lock” on page 53).

NOTICE

You need the right code to do this transaction. You can find the code under section “Code lock” on page 53.

Accessing the OEM menu.

A press on the menu-button takes you to the next menu prompt. After system control has been unlocked, simultaneous presses on the Interval-button and the Nonstop-button call the OEM menu. Following this, the "OEM menu Press Enter" prompt is displayed. When this is confirmed with Enter, the first menu prompt can be worked on. To simplify changing settings, press the UV-button to select the individual number in the numerical value which you want to change. Now use the arrow buttons to enter the wanted number from 0 to 9 at that selected position. Press Enter to save changes only.

The display shows:



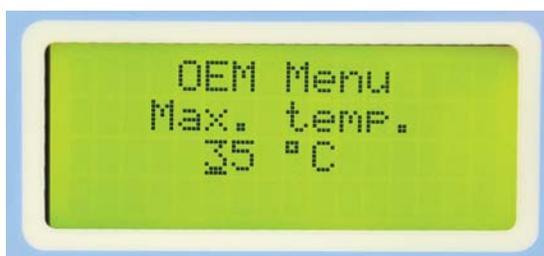
Set the limiting value for temperature

The maximum operating temperature limit for the system is set here. Should this temperature be exceeded, the fault message “Max. temperature” is triggered. This is shown in the 3rd line of the display.

Basic setting: 35 °C

Setting range: 1 - 50 °C

After enter the OEM menu press the menu-button once then the display shows:



Set the rinsing time:

In this Menu point you can set the manually rinsing time.

Basic setting: 30 sec.

Setting range: 10 - 60 sec.

After entering the OEM menu press the menu-button twice then the display shows:



Change the disinfection time

The disinfection time can be set in this menu.

Basic setting: 30 min.

Setting range: 15- 90 min.

After entering the OEM menu press the menu-button 3 times then the display shows:



Set the interval pump time

The interval pump time is the amount of time the pump is working to recirculate water in the system. The standard setting is 5 minutes of pump recirculation for every 30 minutes that the system stands still during Interval mode. The majority of systems do not need this setting to be changed.

Basic setting: 5 min.

Setting range: 1- 30 min.

After entering the OEM menu press the menu-button 4 times then the display shows:



Set the interval rinse time

In this point you can set the rinse interval time. When the system operates in the Interval mode the system is going to be rinsing the hoses for 0.5 sec every 30 min.

Basic setting: 0.5 sec.

Setting range: 0.1- 2 sec.

After entering the OEM menu press the menu-button 5 times then the display shows:



Set the real-time clock

The real time clock can be set in this menu.

Basic setting: The actual date

Setting range: 1-12 Month, 1-31 Day, 0-24 h, 0-60 min.

After entering the OEM menu press the menu-button 6 times then the display shows:



Set the sending interval

In this menu the sending interval between transmissions of measured values and error messages to the RS 232 is set.

Basic setting: 1 hours

Setting range: 0.5-12 hours

After entering the OEM menu press the menu button 7 times then the display shows:

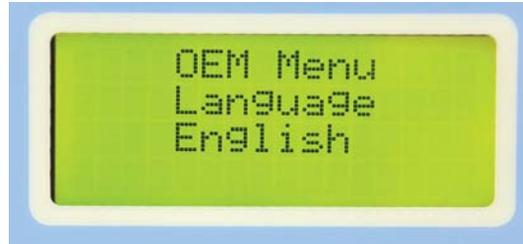


Language selection

Basic setting: English

Setting range: German, English, French

After entering the OEM menu press the menu-button 8 times then the display shows:



Switch units, conductivity/resistance:

Basic setting: Resistance $M\Omega \cdot cm$

Setting range: Resistance $M\Omega \cdot cm$, specific electrical resistance $M\Omega \cdot cm$

After entering the OEM menu press the menu-button 9 times then the display shows:



Switch temperature compensation on/off

Basic setting: On

Setting range: On, Off

After entering the OEM menu press the menu-button 10 times then the display shows:



Printer output

Various parameters are documented by the printer. It differentiates between three types of message:

- Standard message
- Code message
- error message

Standard message:

Here in dependence of the transmit interval of all measured values are printed out. Within the NONSTOP-operation a complete data record is printed out.

Print-out:

e.g.:

```
01.10.10 10:38
GenPure Standard
S.No. 9876/10
Interv. TC on UV off
LF1= 18.2 MΩxcm
LF2= 10.0 MΩxcm
LF3= 0.000 MΩxcm
Temp.= 16.8 °C
TOC= 0 ppb
UV Intens.= 0%
```

The standard record documents all measured values. With systems without TOC measurement and UV-intensity, 0 is entered in place of measured values for these functions!

Code message:

Whenever a code number is entered in system control and confirmed with Enter, the code input is immediately printed out.

Code identification (see “Code lock” on page 62).

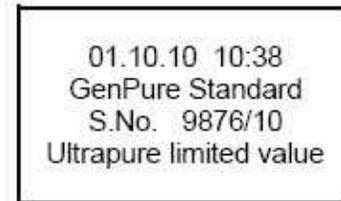
Print-out:

```
01.10.10 10:38
GenPure Standard
S.No. 9876/10
Code 0002
```

Error message:

When a error message is shown in the display, e.g. for the ultrapure water limiting value, then the error message is printed out on expiry of the sending interval.

Print-out:



```
01.10.10 10:38  
GenPure Standard  
S.No. 9876/10  
Ultrapure limited value
```


Maintenance

Contents

- “Maintenance intervals” on page 62
- “Change the ultrapure cartridge” on page 63
- “Disinfection” on page 65
- “Change the ultrafilter” on page 68
- “Structure of the UV-lamp” on page 69
- “Change the UV-lamp” on page 71
- “Change and autoclave the Final filter” on page 75

Regular servicing of your system ensures that the quality of water is maintained. We recommend a service contract with a factory authorized service company to ensure that the system is properly maintained. You then have the certainty of a high operational, safe, and reliable water purification system.

To ensure error-free operation, your system must be checked, serviced and cared for at regular time intervals in accordance with these operating instructions. For this reason, the operating instructions must be readily available to operating and maintenance staff at all times, and be carefully followed.

Calibration of the conductivity is only to be carried out and recorded by a factory-authorized service technician.

Cleaning and disinfection should be performed at least once yearly, or when the ultrapure cartridge is replaced, or when bacteria is present in the product water.



Control and maintenance work on electrical systems are only to be carried out by an appropriately trained, skilled electrician.



Unplug the system from the power outlet for all Maintenance work on the system.

Maintenance intervals

Consumable materials are to be replaced according to the directions below. The intervals were determined for the average user and are completely dependent on the actual feed water quality and volume of water used daily.

Material	Flow chart no.	Catalog no.	Interval	Other problems
Ultrapure cartridge	F1	09.2005	up to 12 Months	Or when the ultrapure water limiting value is exceeded, whichever is shorter. Longer usage can result in bacterial growth on the resin.
Sterile 0.2 micron filter	F2	09.1003	up to 12 Months	Or flow rate is noticeably slower.
Ultrafiltration membrane (only applicable for systems with a UF filter)	F3	50133980	up to 24 Months	Or if there is endotoxin breakthrough in product water or when the water flow rate is markedly slower.
UV-lamp (only applicable for systems with a UV lamp)	UV1	09.2002	up to 24 Months	Or unless system indicates the lamp needs to be replaced.

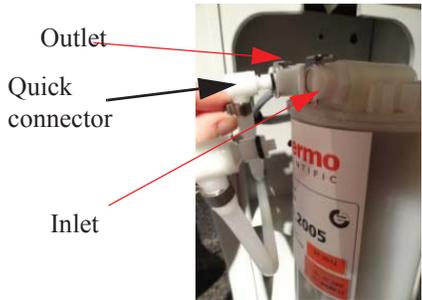
*Please keep in mind that the life of your consumable is directly dependent on the quality of the feed water and the amount of water used daily.

Change the ultrapure cartridge



Replace the ultrapure cartridge when the maximum limiting value that you have set for the ultrapure water is exceeded or when the “change cartridge” message is shown in the display.

Step	Action	Figure
1	Switch the system off and shut off the supply of feedwater.	<p style="text-align: right;">Feedwater supply</p>
2	Remove the cartridge cover by pressing the push button.	
3	Disconnect the quick connectors on the feedwater inlet and purified water outlet of the cartridge and remove the cartridge from the system.	
4	If you change an existing ultrapure cartridge please sanitize your system.	<div style="background-color: #0056b3; color: white; padding: 5px; text-align: center; font-weight: bold; font-size: 1.2em;">NOTICE</div> <p>For sanitize your system see under chapter “Disinfection” on page 65.</p>
5	Locate the new ultrapure cartridge and insert it into the system.	

Step	Action	Figure
6	Plug the quick connects correctly onto the new ultrapure cartridge. You will know they are attached when an audible “click” is heard. Replace the cartridge cover.	

7	Open the supply of feedwater and switch the system on again.	   <p>Feedwater Supply</p>
---	--------------------------------------------------------------	-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

8	<p>NOTICE</p> <p>The Code to do this transaction please refer from the Code table „code lock” found in chapter “Code lock” on page 68. You need a level one code.</p> <ol style="list-style-type: none"> Go in the Menu to the point “change ultrapure cartridge” and press enter. Enter new serial number of the ultrapure cartridge in by pushing the button nonstop or Interval to change the digits and the UV button to go to the next value. When you are finished, press enter and the new serial number is saved. You can only use a serial number one time. 	<p>a. </p> <p>b. </p> <p>c. </p>
---	--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------	----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

9	<p>NOTICE</p> <p>Discard at least 5 liters of water.</p>
---	-----------------------------------------------------------------

Disinfection

CAUTION

Disinfection must be regularly carried out, at the latest when the ultrapure cartridge is replaced, or when bacteria is present in the product water.

Do not stop a disinfection process that is in progress. After faulty disinfection, carry out a new disinfection.

If the system was longer time not in operation, it must be going to be a disinfection process.

A Disinfection cartridge (Catalog no. 09.2201) is required for disinfection of the system.

Use cleaning solutions as follows:

Cleaning Solution, 1 syringe, Catalog no. CMX 25.

NOTICE

For effective disinfection the cartridge must be completely filled with distilled or deionized water.



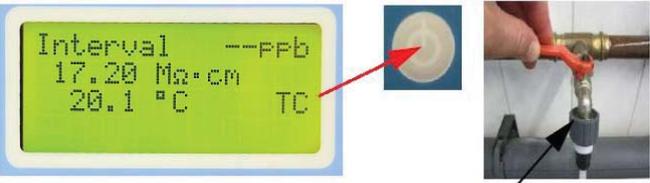
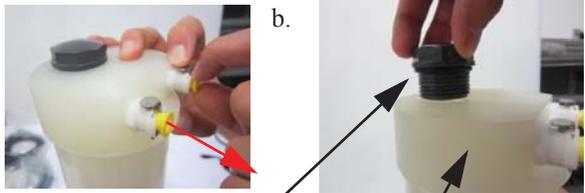
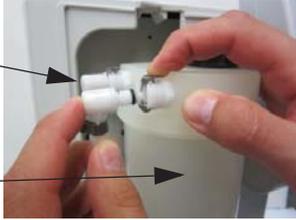
Wear protective gloves for handling syringe of cleaning solution.

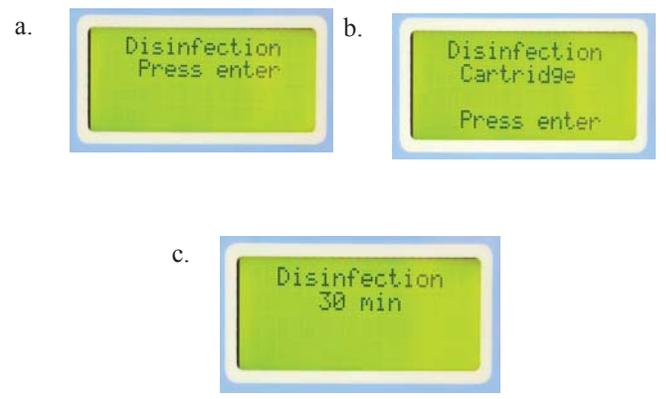
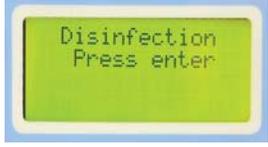
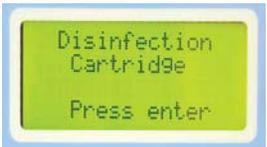
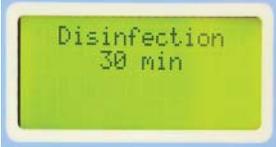


Wear safety goggles when working with cleaning solution.

WARNING

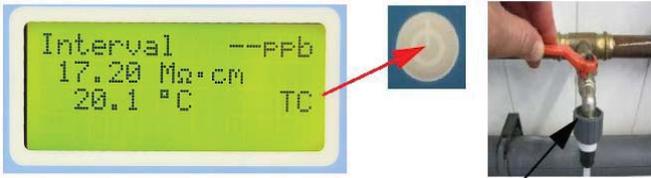
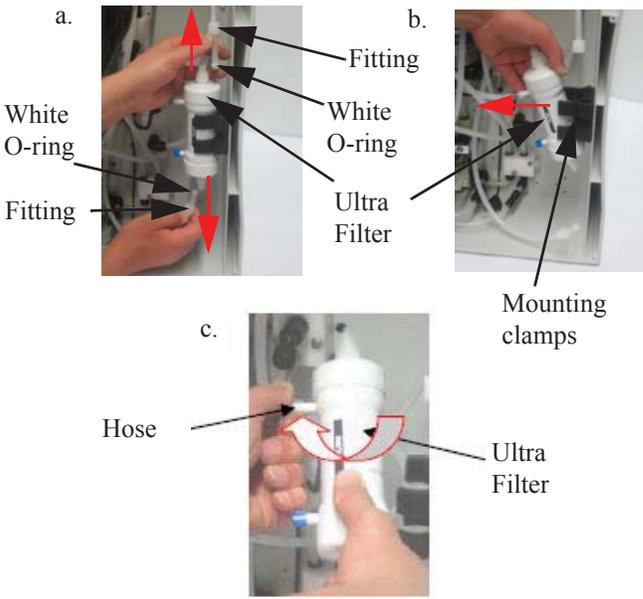
Please observe the information given in the safety data sheet supplied with disinfectant to avoid possible health hazards.

Step	Action	Figure
1	<p>Switch the GenPure System off and shut off the supply of feedwater. After this remove the ultrapure cartridge.</p> <p>NOTICE</p> <p>See under chapter “Change the ultrapure cartridge” on page 63.</p>	 <p>Feedwater supply</p>
2	<p>a. Remove the yellow stoppers.</p> <p>b. Unscrew the stopper from the disinfectant cartridge.</p> <p>c. Fill the cartridge with distilled water then empty the contents of a box respectively a syringe of Cleaning solution into the water.</p>	 <p>Stopper</p> <p>Disinfectant cartridge</p>
3	<p>Screw the stopper back on the disinfectant cartridge and connect the cartridge into the system.</p> <p>NOTICE</p> <p>See under chapter “Change the ultrapure cartridge” on page 63 to put in the ultrapure cartridge in to the system.</p>	 <p>Quick connectors</p> <p>Disinfectant cartridge</p>
4	<p>Re-open the feedwater supply, switch the system on again.</p>	 <p>Feedwater Supply</p>

Step	Action	Figure
5	<p>Push the menu button until “Enter code” is displayed</p> <p>NOTICE</p> <p>The Code to do this transaction please refer from the Code table under chapter “Code lock” on page 68. You need a level 1 code.</p> <ol style="list-style-type: none"> Select “Disinfection” from the system menu and press “Enter”. Confirm the Disinfection Cartridge has been loaded by pushing “Enter” again The disinfection process will begin. <p>NOTICE</p> <p>The disinfection program is finished after approx 30 min and is adjustable in the OEM Menu.</p>	 <p>a. </p> <p>b. </p> <p>c. </p>
6	<p>Switch the system off and shut off the water supply.</p>	<p>NOTICE</p> <p>See step 1.</p>
7	<p>Remove the disinfectant cartridge, empty and dry it and put in the yellow stoppers that you have saved for later use. Save the disinfection cartridge for later use.</p>	<p>NOTICE</p> <p>See step 5 under chapter “Change the ultrapure cartridge” on page 63.</p>
8	<p>CAUTION</p> <p>Before dispensing water from the system, let water run out for approx 15 minutes. The system is then ready for use.</p>	

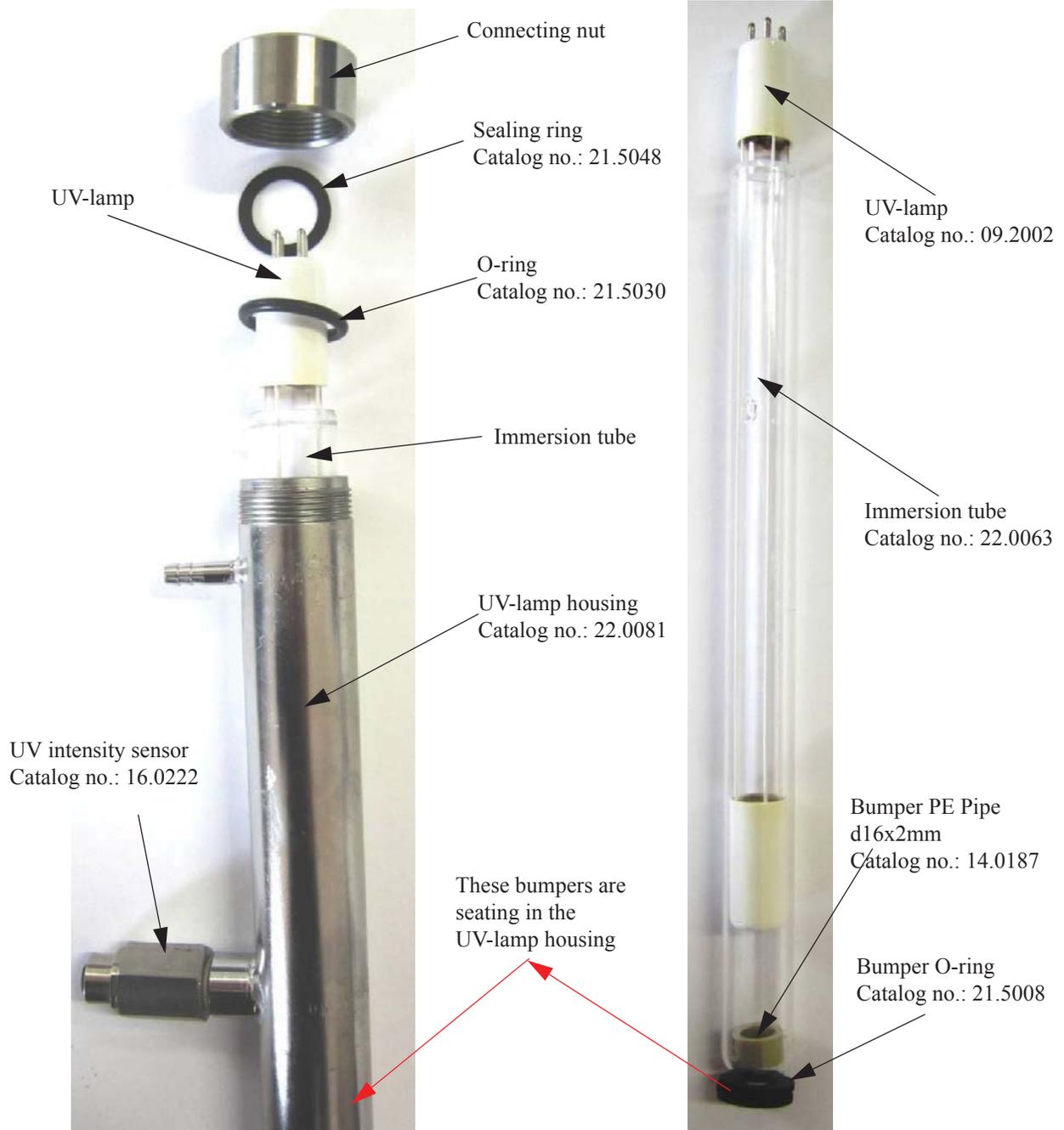
Change the ultrafilter

(applicable only for systems with UF)

Step	Action	Figure
1	Switch the GenPure System off and shut off the supply of feedwater.	 <p>Feedwater supply</p>
2	<p>Remove the four screws of the back panel.</p> <p>CAUTION</p> <p>Remove carefully the back panel from the system and unscrew the yellow ground wire from the back panel.</p>	 <p>Screws</p>
3	<p>a. Pull out the hoses 8 mm by unscrew the two fittings (see red arrows).</p> <p>b. After this procedure draw out the ultrafilter from the mounting clamp (see red arrow).</p> <p>c. Hold with one hand the hose and with the other hand turn in clockwise direction the ultrafilter to unscrew the hose connection.</p> <p>d. When you are finished with step c install the new ultrafilter by attaching hoses and mounting it in the clamp.</p> <p>NOTICE</p> <p>e. When you are installing the new UF filter the flow arrow of the filter must be pointing to the bottom of the system.</p>	 <p>a. Fitting, White O-ring, Fitting</p> <p>b. Fitting, White O-ring, Ultra Filter, Mounting clamps</p> <p>c. Hose, Ultra Filter</p>
4	Reinstall the back panel, reopen the feedwater supply and switch on the system again.	 <p>Feedwater Supply</p>

Structure of the UV-lamp

UV unit with UV intensity sensor



UV unit without UV intensity sensor



Change the UV-lamp

⚠ WARNING

Never look directly into a switched-on UV-lamp, as UV-light endangers eyesight!



Always wear safety gloves when changing the UV-lamp, in order to prevent that your skin comes in contact with the UV-lamp glass.



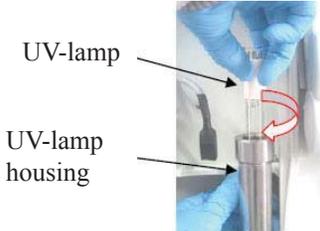
Wear directly a breathing protector when you are seeing that the glass of the UV-lamp is broken and ventilate the room well.

NOTICE

Contact your local Service organization to proceed as the proper disposal of the used UV-lamp. The Hg content in the UV-lamp is so low so that no damage to the environment can arise.

(applicable only for systems with UV lamp)

Step	Action	Figure
1	Switch the GenPure System off and shut off the supply of feedwater.	
2	Remove the cartridge cover and take off the ultrapure cartridge.	See under chapter “Change the ultrapure cartridge” on page 63.
3	Unscrew the bracket from the mounting plate and take it up over the UV-lamp cable.	
4	Draw the UV-lamp housing slightly to the front (see red arrow) to the front and take the plug off of the UV-lamp.	

Step	Action	Figure
5	Now carefully draw the UV-lamp upwards while lightly turning it clockwise. During the replacement of a UV-lamp, great care must be taken to avoid touching the glass of the UV-lamp with fingers, to avoid dirtying of the lamp which would impair the functioning of it.	 <p>UV-lamp</p> <p>UV-lamp housing</p>
NOTICE		NOTICE
We therefore recommend that clean gloves be worn.		See chapter “Structure of the UV-lamp” on page 69 where is seating the sealing ring to not damage it.

Step	Action	Figure
------	--------	--------

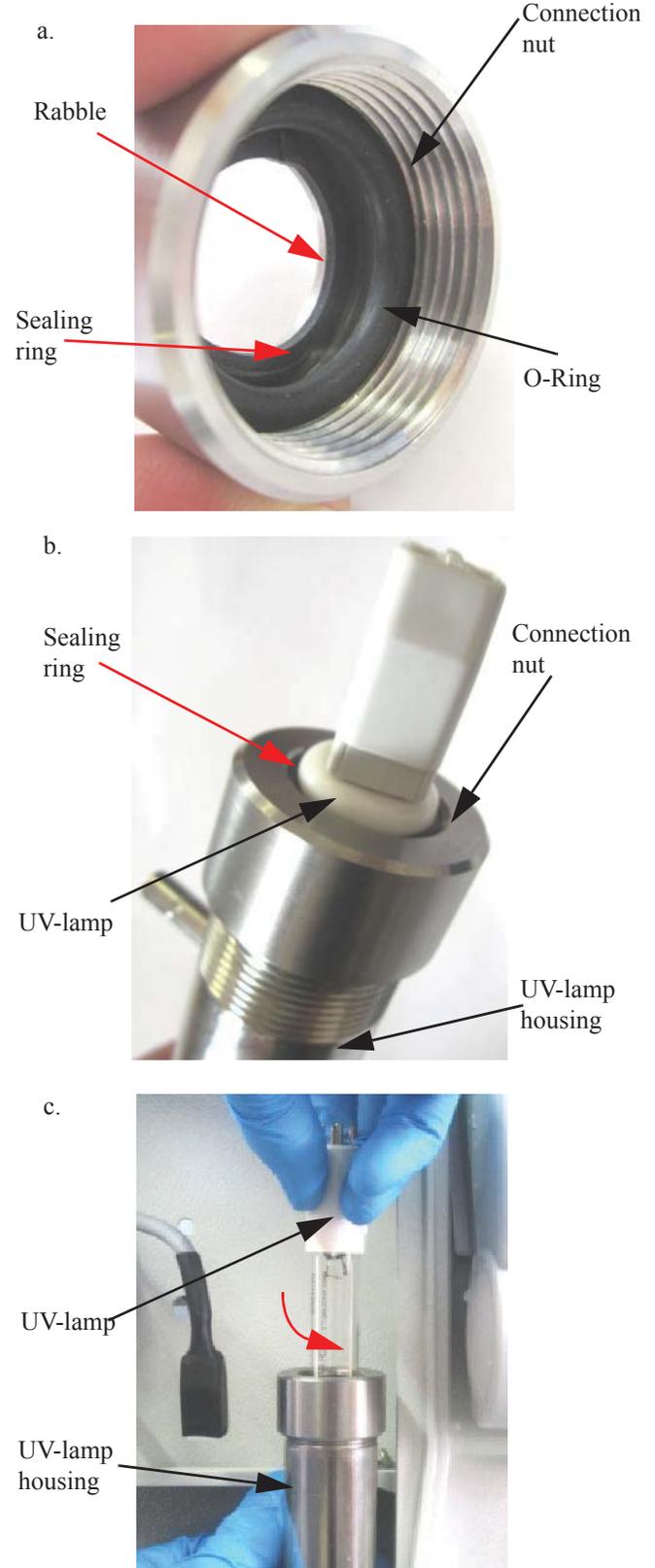
6

CAUTION

Ensure that the position of the sealing ring (flat o-ring at the top of the connection nut) is correct as you put in the new UV-lamp, otherwise you will have a leak. The sealing ring must be seat in the rabbet of the connecting nut (see picture a and b).

Carefully introduce the new UV-lamp under a slight turning motion like before but in the anti-clockwise direction (see picture c). Attach the plug into the lamp and push the housing back to the system. Once it is in place, re-mount the bracket holding the UV housing onto the system's the mounting plate.

Figure



Step	Action	Figure
7	Put the cartridge cover back on (see under section “Change the ultrapure cartridge” on page 63), re-open the feed water supply and switch the system on again.	 Feedwater Supply

8 Push the menu button until “Enter code” is displayed.

NOTICE

The Code to do this transaction please refer from the Code table under section “Code lock” on page 68. You need a level 3 code.

- a. After entering the code push the Menu and UV button simultaneously. The display shows UV Menu.
- b. Push the Menu button repeatedly until new UV-lamp appears and press enter to confirm.
- c. The system sets the operating hours counter of the UV-lamp back and save the new values by an automatically calibration.

NOTICE

The UV-lamp must be switched on (Nonstop mode).
The calibration process of the UV intensity can be take between 5 min. and 2 hours.

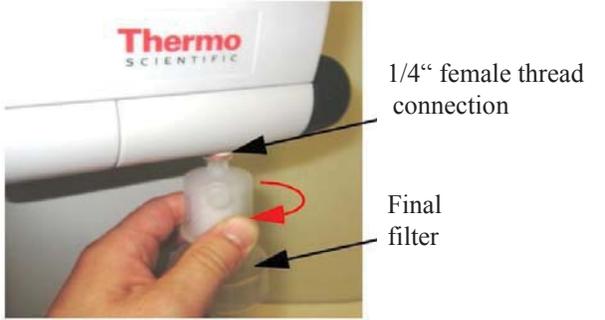
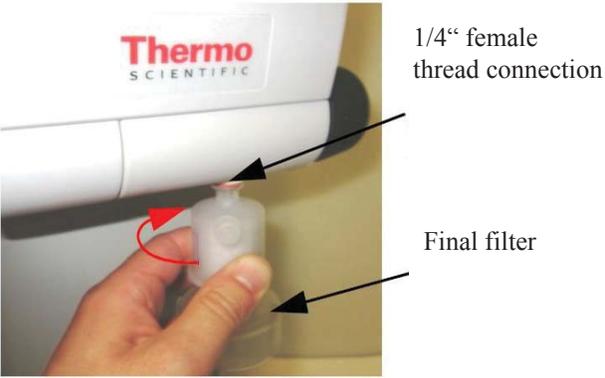
a.



b.



Change and autoclave the Final filter

Step	Action	Figure
1	Screw out the blocked or used final filter by turn it in clockwise direction.	
2	Unpacking the new Final filter and screw in the filter counter clockwise (See arrow in the picture) in the dispensing valve outlet (R 1/4" female thread)	
3	Rinse about 3 liters through final filter before use.	

Autoclave the Final filter

NOTICE

To increase the lifetime of the filter you can autoclave it. To autoclave the final filter proceed as follows.

Step	Action	Figure
1	Unscrew the used final filter by turn it in clockwise direction.	

2 Use a autoclave to sterilize the filter.

The temperature of the autoclaving process must be 121°C and should take 30 min. You can repeat the procedure for the filter up to 5 times. When the sterilization is finished screw in the final filter back to the dispensing valve outlet (see chapter “[Change and autoclave the Final filter](#)” on [page 75](#)).

NOTICE

If you are trying to dispense water and nothing is coming out from the outlet, the final filter is blocked. Please look then in chapter “[Trouble shooting](#)” on [page 79](#) or change with a new one.

Waste disposal

NOTICE

Before you are shipping your Thermo Scientific Barnstead Ultrapure Water System to the waste disposal, only a specially trained personal is to be taken out the system of operation.

When the packaging is no longer needed it can be disposed of as household waste.

Systems are in conformity with EEC Guideline 2011/65/EU.

The system is not to be thrown away as household waste but must be properly disposed of. It can be returned to the manufacturer for safe disposal according to EEC Guideline 2011/65/EU. We therefore request our customers in Germany and other member States in the European Economic Area to contact our local service center or our headquarters or per E-Mail to:

weee.recycle@thermofisher.com

WEEE-Reg.-no.: DE 12471402

In countries outside of the European Economic Area, please contact your local authorities or waste disposal company.

Trouble shooting

NOTICE

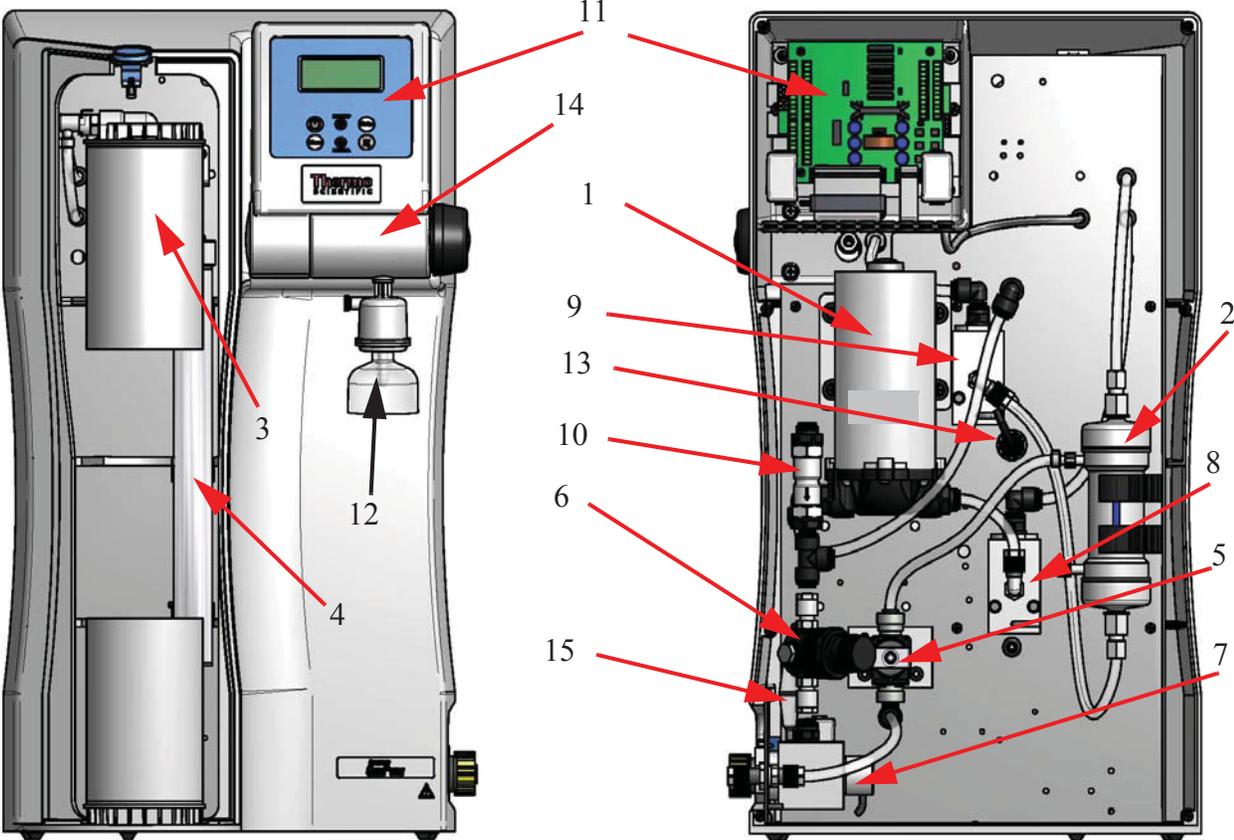
If the error can not be solved by the customer, the service is should be to refrain.

Error	Cause	Remedy
The system does not start	<ul style="list-style-type: none"> No supply of power 	<ul style="list-style-type: none"> Provide power
Dispensing not possible	<ul style="list-style-type: none"> Feedwater tap is closed Feedwater and rinse water connections are mixed up Feedwater pressure < 0.1 bar Final Filter is blocked 	<ul style="list-style-type: none"> Open the feedwater tap Correct the connections Increase the feedwater pressure Change with a new one
Resistance < 18.2 MΩ·cm	<ul style="list-style-type: none"> Ion exchange capacity is exhausted Poor feedwater Temperature compensation turned off calibration needed 	<ul style="list-style-type: none"> Replace cartridge with a new one Correct feedwater Turn temperature compensation on (Display should show “TC” in bottom right) Contact Service for calibration
System control no longer reacts	<ul style="list-style-type: none"> Improper operation error PCB Faulty Dispense button 	<ul style="list-style-type: none"> Unplug the mains plug for 5 seconds. Contact the Service. Contact Thermo for service
Water flows out	<ul style="list-style-type: none"> Leaky hose connection Feedwater pressure > 6 bar 	<ul style="list-style-type: none"> Check and seal the hose connection Install a pressure reducer Contact Thermo for service

Error	Cause	Remedy
Dispensed amount is too small	<ul style="list-style-type: none"> • UF-Module blocked • Feed water pressure too low • Internal pressure too low • Volumetric Dispense out of Tolerance 	<ul style="list-style-type: none"> • Replace UF-module • Increase the feed water pressure • Readjust pressure reducer • Contact Thermo for volume calibration
Wrong time or date	<ul style="list-style-type: none"> • Time zone • Summer/winter time 	<ul style="list-style-type: none"> • Reset time and date
Wrong language	<ul style="list-style-type: none"> • Wrong language set 	<ul style="list-style-type: none"> • Correct the language setting
Error message: <i>"Limit value feed"</i>	<ul style="list-style-type: none"> • Feedwater conductivity too high • Limiting value set too low • TOC selected on non-TOC units 	<ul style="list-style-type: none"> • Check the pretreatment • Check and suit the limiting value setting • Turn LF3 to off
Display reads +IN	<ul style="list-style-type: none"> • Measuring cell cable break 	<ul style="list-style-type: none"> • Replace measuring cell
Error message: <i>"Lim. va.pure w."</i>	<ul style="list-style-type: none"> • Ultrapure cartridge exhausted • Limiting value set too low 	<ul style="list-style-type: none"> • Replace with new ultrapure cartridge • Check and set the limiting value
Error message: <i>"UV-duration"</i>	<ul style="list-style-type: none"> • UV-Lamp operating time has been exceeded 	<ul style="list-style-type: none"> • Replace the UV-lamp • Re-set the operating time counter
Error message: <i>"UV-intensity"</i>	<ul style="list-style-type: none"> • UV-Lamp intensity no longer sufficient • UV-Sensor is dirty • Limiting value set too low 	<ul style="list-style-type: none"> • Replace and measure in a new UV-lamp • Clean the UV-sensor • Check and set the limiting value
Error message: <i>"max. Temperature"</i>	<ul style="list-style-type: none"> • The temperature in the system is too high • Interval pump time too long • Limiting value set too low • Feedwater temperature is too high 	<ul style="list-style-type: none"> • Reduce the temperature by running water off • Reduce interval pump time • Check and suit the limiting value • Reduce the feedwater temperature
Error message: <i>"Measuring cell LF1"</i>	<ul style="list-style-type: none"> • Measuring cell cable break • System control defect • Conductivity of ultrapure water outside of the measuring range 	<ul style="list-style-type: none"> • Replace the measuring cell • Replace system control • see "Resistance < 18.2 MW·cm" on page 79

Error	Cause	Remedy
Error message: "Measuring cell LF2"	<ul style="list-style-type: none"> • Measuring cell cable break • System control defect • Feedwater conductivity outside of measuring range 	<ul style="list-style-type: none"> • Replace the measuring cell • Replace system control • see "Error message: "Limit value feed"" on page 80
Error message: "Measuring cell LF3"	<ul style="list-style-type: none"> • Measuring cell cable break • System control defect 	<ul style="list-style-type: none"> • Replace the measuring cell • Replace system control
Error message: "Temp. meas. cell."	<ul style="list-style-type: none"> • A break in the measuring cell cable • System control defect 	<ul style="list-style-type: none"> • Replace the measuring cell • Replace the system control
Error message: "change cartridge"	<ul style="list-style-type: none"> • Operating hours of the filter cartridge has expired 	<ul style="list-style-type: none"> • Replace it with a new one

Replacement parts



Parts marked with an „x“ must be changed by a qualified service professional

Pos.	Designation	Catalog no.	
1	Pressure boosting pump	50149262	x
2	Ultrafiltration module (optional)	50133980	
3	Ultrapure cartridge	09.2005	
4	Replacement UV-lamp	09.2002	
5	Rinsing solenoid valve	50131190	x
6	Pressure reducer	50133985	
7	Feedwater conductivity measuring cell	16.0126	
8	Ultrapure water conductivity measuring cell	50133992	
9	TOC conductivity measuring cell (optional)	50133992	
10	Check valve	50150598	
11	Microprocessor system control CPU board	50131348 26.0025	
12	Final filter 0.2µm, for dispensed water	09.1003	
13	UV-Intensity sensor (optional)	16.0222	
14	Dispensing valve	50133988	x
15	Fuse holder for glass tube fuse 5 x 20 mm Glass tube fuse 5 x 20 mm, 3,15 A, slow fuse	50143154 50150714	
16	Tabletop power pack, 24V DC (not shown)	50149597	
17	UV- ballast (not shown)	22.0088	

NOTICE

We ask for your understanding that our guarantee for this system is invalidated when replacement parts, accessories or consumable materials from other manufacturers are used in or for the system, as we have no influence on their composition or quality.

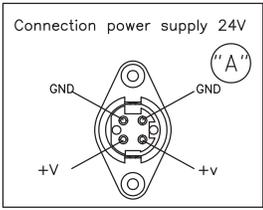
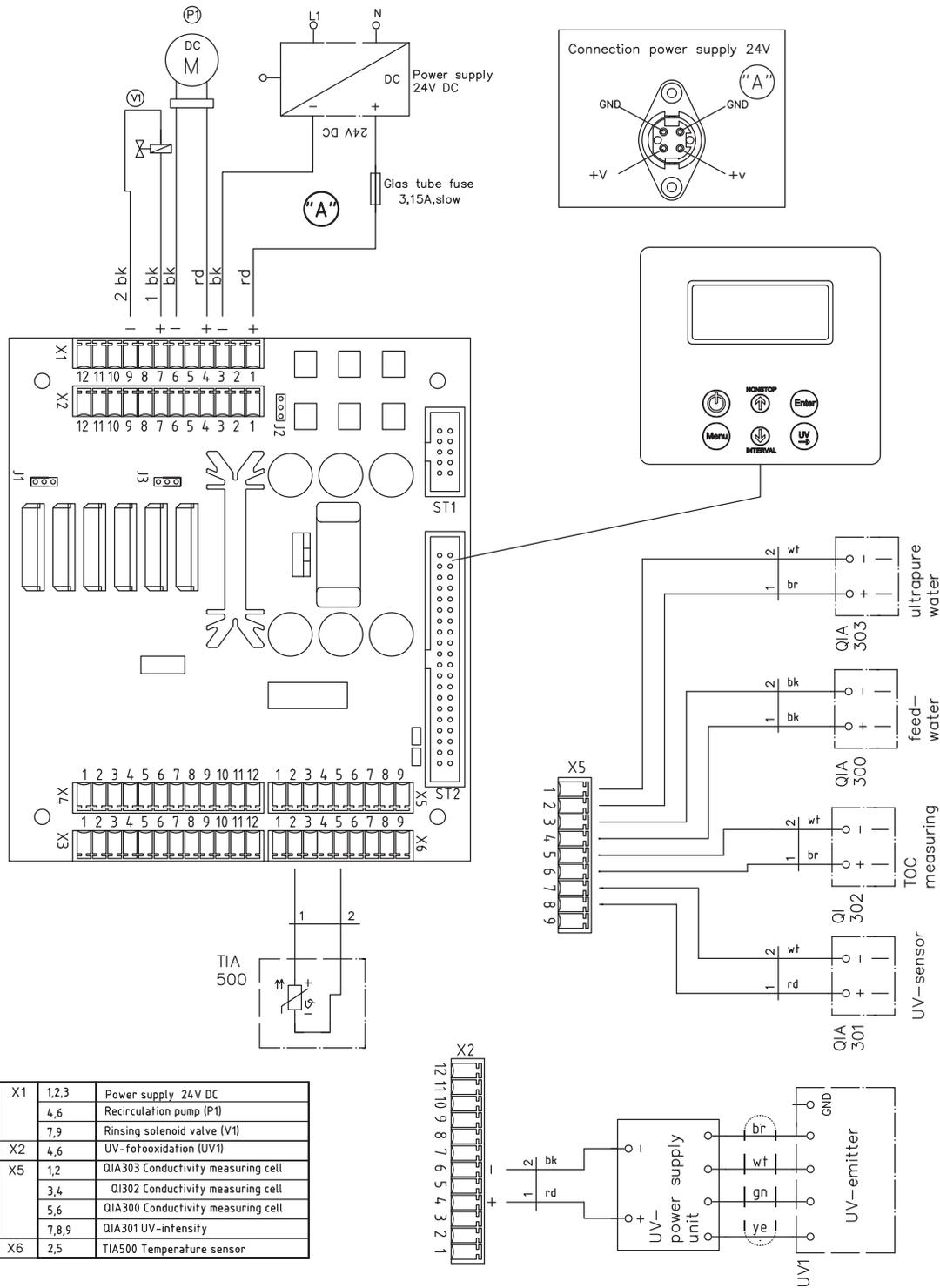
Consumable materials

Designation	Catalog no.
Ultrapure cartridge	09.2005
UV-Lamp	09.2002
Ultrafiltration module	50133980
Final filter 0.2 μm	09.1003

Accessories

Designation	Catalog no.
Disinfection cartridge	09.2201
Cleaning Solution, 1 syringe	CMX25
Printer	09.2207
Ion exchanger DI 1500	02.1500
DI 1500 hose kit for new installations	04.1690

Terminal assignments



X1	1,2,3	Power supply 24V DC
	4,6	Recirculation pump (P1)
	7,9	Rinsing solenoid valve (V1)
X2	4,6	UV-fotooxidation (UV1)
X5	1,2	QIA303 Conductivity measuring cell
	3,4	QI302 Conductivity measuring cell
	5,6	QIA300 Conductivity measuring cell
	7,8,9	QIA301 UV-intensity
X6	2,5	TIA500 Temperature sensor

Maintenance records

Customer address: _____ **Location:** _____

_____ **System type:** _____

_____ **Serial no.:** _____

_____ **Year made:** _____

Date	Feedwater resistance [MΩxcm]	Ultrapure water resistance [MΩxcm]	Temperature [°C]	TOC value [ppb]	UV intensity [%]	UV-lamp operating time [h]

Ultrapure water flow rate [l/h]	Last filter cartridge replacement	Last cleaning, disinfection	Remarks	Signature

Any false entry is considered to be a falsification of documents.

The following point should be observed for maintenance of the quality of the system:

- 1x / Weekly, acquire measured values.

Contact Information Thermo Scientific

The address to contact when your system requires service:

Overview of Thermo Scientific International Sales Organization

Postal address USA:

Thermo Scientific
275 Aiken Road
Asheville, NC 28804
USA

Enquiries from USA/Canada

Sales: +1 866 984 3766

Service: +1 800 438 4851

Enquiries from Latin America

Sales: +1 866 984 3766

Service: +1 866 984 3766

Enquiries from Asia:

China

Sales: +86 10 8419 3588

Service : Toll free 8008105118
Support Mobile 4006505118 or +86 10 8419 3588

India

Sales: +91 22 6716 2200

Service: Toll free 1 800 22 8374 or +91 22 6716 2200

Japan

Sales: +81 45 453 9220

Service: +81 45 453 9224

Enquiries from the Rest of Asia/Australia/New Zealand

Sales: +852 2885 4613

Service: +65 6872 9720

Enquiries from Countries not listed / Rest of EMEA

Sales: +49 6184 90 6940 or +33 2 2803 2000

Service: +49 6184 90 6940

Enquiries from Europe:

Austria

Sales: +43 1 801 40 0

Service : +43 1 801 40 0

Belgium

Sales: +32 53 73 4241

Service: +32 53 73 4241

Finland/Nordic/Baltic countries

Sales: +358 9 329 100

Service: +358 9 329 100

France

Sales: +33 2 2803 2180

Service : +33 825 800 119

Germany:

Postal Address Germany:

Thermo Electron LED GmbH

Robert-Bosch-Straße 1

D - 63505 Langenselbold

Phone

Sales Toll free 0800 1 536 376

or +49 6184 90 6940

Service Toll free 0800 1 112110

or +49 6184 90 6940

E-Mail info.labequipment.de@thermofisher.com

Italy

Sales +39 02 95059 341

Service +39 02 95059 250

Netherlands

Sales +31 76 579 5555

Service +31 76 579 5639

Russia/CIS

Sales +7 812 703 4215

Service +7 812 703 4215

Spain/Portugal

Sales +34 93 223 0918

Service +34 93 223 0918

Switzerland +41 44 454 1212

Service +41 44 454 1212

UK/Ireland

Service +44 870 609 9203

Sales +44 870 609 9203

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

Index

A

- Accessories 87
- Airborne sound emission 19
- Ambient conditions 19
- Assembly of the holder of the dispenser 27
- Autoclave
 - Final filter 76
- Autoclave the Final filter 76
- Available GenPure Pro Systems 14

C

- Cell constants of the measuring cells 19
- Change
 - Final filter 75
 - Ultrafilter 68
 - UV-lamp 71
- Change and autoclave the Final filter 75
- Change the ultrapure cartridge 63
- Change the UV-lam 71
- Code message 58
- Complaints 6
- Connectors for water 19
- Consumable materials 85
- Contact information 93
- Contact Information Thermo Scientific 93
- Contents 1

D

- Description of Display 40
- Dimension and weight 18
- Disinfection 51, 65

E

- Electrical connections / external switched mode power supply 19
- Entering a code number 53
- Error history 52
- Error message 59
- Examination on receipt 6
- Explanatory notes v
- Explanatory notes on the operating instructions v
- Extent of assembly kit 12
- Extent of delivery 11

F

- Fault message 59
- Filter cartridge
 - Change 63
- Final
 - Autoclave 76
- Final filter
 - Change 75
- Flow chart control unit 40
- Flow chart of menu system control 41
- Flow chart 33

H

- Holes for wall mounting 28
- How the system functions 35

I

- Information, legal iv
- Installation 23
- Installation area 21
- Installation examples 31
- Intended use 15
- Interval operating mode after switching on 44
- Interval operation 45

L

- Legal iv
- Legal Information iv

M

- Maintenance 61
 - Intervals 62
 - Records 91
- Maintenance intervals 62
- Maintenance records 91
- Materials of parts which contact water 20
- Mounting the power pack 29

N

- Non-stop mode 45

O

- OFF mode 46

Index:

- Operating elements 39
 - Operating modes 44
 - Interval operating mode after switching on 44
 - Interval operation 45
 - Non-stop mode 45
 - OFF mode 46
 - UV-Lamp 45
 - P**
 - Packaging 5
 - Packing for return shipment 6
 - Preface iii
 - Print out of Data 52
 - Printer output 58
 - Print-out 58
 - Product water quality 17
 - Putting into operation 37
 - Putting system into operation 37
 - R**
 - Replacement parts 83
 - Return shipment 6
 - S**
 - Safety precautions 7
 - Specifications 17
 - Standard message 58
 - Standards and Directives vi
 - Start your system into operation 25
 - Structure of the UV-lamp 69
 - System
 - Functions 35
 - Putting into operation 37
 - System control 43
 - General information 44
 - T**
 - Technical 17
 - Technical specifications 17
 - Airborne sound emission 19
 - Ambient conditions (DIN EN 61010-1 (VDE 0411-1) 2011-02) 19
 - Cell constants of the measuring cells 19
 - Connectors for water 19
 - Demands the feedwater must fulfill 17
 - Dimension and weight 18
 - Electrical connections / external switched mode power supply 19
 - Materials of parts which contact water 20
 - Product water quality 17
 - Terminal assignment 89
 - Terminal assignments 89
 - The installation area 21
 - Transport 5
 - Transport and packaging 5
 - Trouble shooting 79
- U**
 - Ultrafilter
 - Change 68
 - Ultrapure water limiting value 48
 - Unintended Use 15
 - User menu 47
 - Disinfection 51
 - Entering a code number 53
 - Error history 52
 - Feedwater measured value and limiting value 47
 - Print out of Data 52
 - Registering the station 53
 - Rinsing the ultrafilter 50
 - Ultrapure water limiting value 48
 - UV Intensity and operating time 48
 - UV-Lamp 46
 - UV-lamp
 - Change 71
 - Structure 69
 - UV-Lamp operating time and intensity 49
 - W**
 - Wall mounting
 - Holes 27
 - Warranty iv
 - Waste disposal 77

Index:



thermoscientific.co

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. 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
81 Wyman Street
Waltham, MA 02451

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific