



Thermo Fisher Scientific Barnstead GenPure with xCAD Ultrapure water system

Operating instruction

50131281 Revision A November 2013

© 2013 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

© 2013 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

NOTE

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 GenPure with xCAD ultrapure Water System 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 and accessories that have been approved by Thermo Electron LED GmbH are used (third-party spares 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.



Important operating and/or maintenance instructions. Read the operating instructions with due care.

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.

The information provided in these operating instructions is only valid for the system which has the serial number which is to be entered on the front page. This information is valid for the system that is received.

NOTE

Please enter the serial number* of your GenPure with xCAD system in the space provided on the front page.

* Read the serial number of your ultrapure water system from the type plate.

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
- The catalog number

Standards and Directives

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

- Low Voltage Directive 2006/95/EC
- EMC Directive 2004/108/EC
- Machine directive 2006/42/EC
- ASTM D1193-6

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 Information	iv
Warranty	iv
Explanatory notes on the operating instructions	v
Standards and Directives	vi
 Chapter 1 Transport and packaging	5
Examination on receipt	6
Complaints	6
Packaging and return shipment	6
 Chapter 2 Safety precautions	7
 Chapter 3 Extent of delivery	9
Extent of assembly kit	10
Available GenPure with xCAD versions	12
GenPure with xCAD bench version:	12
GenPure with xCAD wall version:	13
 Chapter 4 Intended use	15
 Chapter 5 Technical specifications	17
 Chapter 6 The installation area	23
 Chapter 7 Installation	25
Connectors GenPure system	26
Connectors xCAD	27
Installation of GenPure with xCAD system, bench version	28
Installation of GenPure with xCAD system, wall version	32
Wall mounting GenPure system	37
Mounting the power pack (voltage supply)	39
Installation examples	41
 Chapter 8 Flow charts	42
Flow chart, GenPure Standard with xCAD	43
Flow chart, GenPure UV with xCAD	44
Flow chart for GenPure UF	45

Flow chart, GenPure UV/UF with xCAD	46
Flow chart, GenPure UV-TOC with xCAD	47
Flow chart, GenPure UV-TOC/UF with xCAD	48
Chapter 9 How the system functions	50
Chapter 10 Putting system into operation	52
Chapter 11 Operating elements.....	54
Flow chart control unit	55
Flow chart control unit	56
Chapter 12 The system control	58
General information	59
Operating modes	59
Interval operating mode after switching on	59
Non-stop mode	60
Interval operation	60
UV lamp	60
Water dispensing via volumetric dispense	61
OFF mode	62
User menu	62
Feedwater limiting value:	62
Ultrapure water limiting value	63
UV-Lamp operating time and intensity:	63
Ultrapure cartridge operating hours counter:	64
Rinsing procedure	64
Disinfection procedure	65
Error history	66
Print out of Data	66
Entering a code number	67
Code lock	67
The OEM Menu	68
Set the limiting value for temperature:	68
Set the rinsing time:	68
Change the disinfection time:	69
Set the pump interval time:	69
Set the rinse interval time:	70
Adjust the real time clock:	70
Set the sending interval:	70
Select the language:	71
Switch units, conductivity/resistance:	71
Switch temperature compensation on/off:	71
Using volume control for water dispensing	72
Printer output	72
Standard message:	72
Code message:	73
Error message:	73

Chapter 13	Maintenance	74
	Maintenance intervals	75
	Change the ultrapure cartridge	76
	Disinfection	78
	Change the ultrafilter	81
	Structure of the UV lamp	83
	Change the UV lamp	85
	Change and autoclave the Final filter	89
	Autoclave the Final filter	90
Chapter 14	Waste disposal	92
Chapter 15	Trouble shooting	94
Chapter 16	Replacement parts and consumables	98
	GenPure	98
	xCAD	100
Chapter 17	Consumable materials	102
Chapter 18	Accessories	104
Chapter 19	Terminal assignment	106
Chapter 20	Maintenance records	108
Chapter 21	Contact Information Thermo Scientific	110

Transport and packaging

Contents

- “Examination on receipt” on page 6
- “Complaints” on page 6
- “Packaging and return shipment” on page 6

Ultrapure water systems are carefully checked and packed prior to shipment but, despite this, there is always a possibility that damage could occur to them during shipment

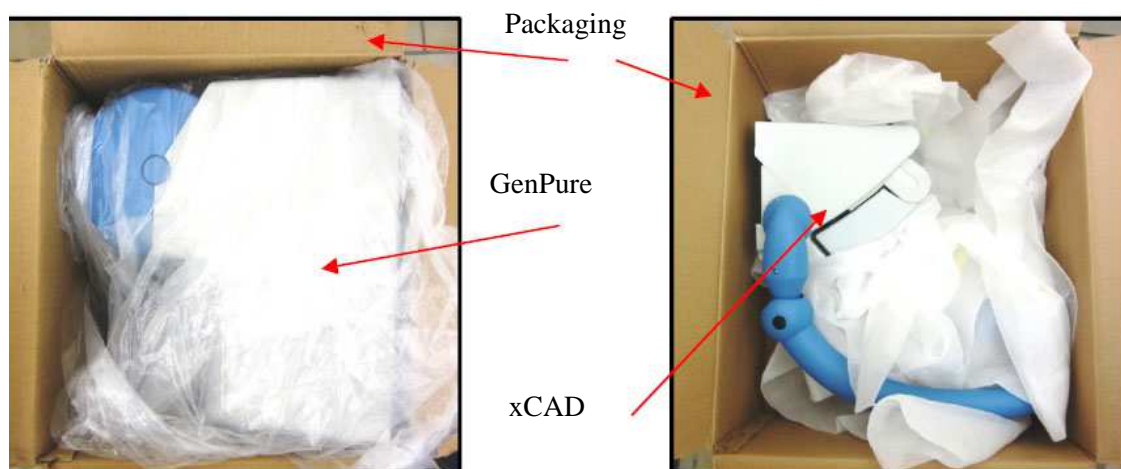
Examination on receipt

- Check the completeness of the delivery against the shipping papers.

NOTE

Does the packaging show signs of damage?

- Inspect the equipment supplied for damage.



Complaints

Should the system have been damaged during transport:

- Immediately contact your delivery transport agency.
- Keep the packaging, including the outer cardboard box, for a possible inspection and/or return shipment.

Packaging and return shipment

Whenever possible use the original packaging and packing materials.

Should these no longer be available:

- Pack the system in packing film and then in a strong cardboard box so that it is held shock-proof.

NOTE

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.

Safety precautions

NOTE

For your own safety, please observe the above safety precautions.



DANGER

Your GenPure series system is a contemporary ultrapure water system. It serves exclusively to purify pretreated water. The water it produces is not fit for drinking.

- 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.
- Note that the manufacturer is not liable for damages that result from improper operation of the system, or from use of it for other than the intended purpose.
- 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.
- Protective means need to be installed to prevent tap water contamination.
- A grounded 100-250V, 50/60Hz socket must be available.
- The installation area must have a drain at floor level with at least a nominal diameter of 63 mm or 2.48 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.
- 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 75.
- The surface or wall where the system is to be installed must have sufficient load-bearing capacity (see „Technical specification“ on page 18).
- When installing the ultrapure water system, ensure that there is sufficient working area around it for convenient operation of it (e.g. ultrapure cartridge replacement, connections, etc.)...



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



To avoid possible risks of crush injury, cuts or electric shock when handling the system, never take the protective casing off of the system. Only trained, skilled personnel are to be assigned to carry out maintenance of the system.

- Regularly carry out visual inspection of the system before operating it, as splashes of liquid could result in a danger of slipping. Any emergent liquid must be immediately mopped up.



Wear protective gloves when chlorine tablets or a disinfection syringe (only US) are to be handled during maintenance. Do not stop a disinfection process that is in progress. After faulty disinfection, carry out a new disinfection run.



Increased heat might be caused by system or system component defects. To reduce skin damage it is recommended to wear protective gloves.

- Do not use oxidative cleaning agents when cleaning the system. They would cause damage to it.
- If the system has a defect, proceed as follows:
 - Switch the system off (dead)
 - Stop the water inlet
 - Contact the Local service organization

Extent of delivery

Contents

- “Extent of assembly kit” on page 10
- “Available GenPure with xCAD versions” on page 12

Extent of assembly kit

Ultrapure cartridge
Catalog no.: 09.2005



Final filter 0.2 µm
Catalog no.: 09.1003

NOTE

To increase the lifetime of the filter a sterilization at 120 °C for 30 min is recommended. The procedure for the filter can be repeated up to 10 times.



Transformer-table power pack
Catalog no.: 50134196



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



Feedwater connecting kit
Catalog no.: 25.0075



PE hose, Ø8mm x 2 m or 0.31 inch x 2.19 yard
Catalog no.: 18.0036



PE hose, Ø6mm x 5 m or 0.24 inch x 5.46 yard
Catalog no.: 18.0047



PE hose, Ø4mm x 5 m or 0.16 inch x 5.46 yard
Catalog no.: 18.0053



Connecting Cord (US) Catalog no.: 50132200
Connecting Cord (british) Catalog no.: 50132203
Connecting Cord (euro) Catalog no.: 50132215



Mounting parts for wall mounting GenPure system and xCAD
wall version:

-Plug 4 x S6

Catalog no.: 21.0002 (for xCAD)

-Screw 4x40 mm or 4 x 1.57 inch

Catalog no.: 21.0001 (for xCAD)

-Plug 2 x S8

Catalog no.: 21.0035 (for GenPure)

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

Catalog No.: 21.0057 (for GenPure)



3 Extent of delivery

Available GenPure with xCAD versions

2x Sub-D extension cable,
25 pin, 5 m or 5.46 yard
Catalog no.: 16.0375



Spiral cable, inlet Ø 9 mm or 0.35 inch,
B10 – 40 mm or B10 - 1.57 inch
Catalog no.: 16.0340



Available GenPure with xCAD versions

GenPure with xCAD bench version:



50131286 standard
50131250 UF
50131254 UV
50131252 UV/UF

50131296 UV-TOC

50131298e UV-TOC/UF

Basic system
Basic system + ultrafiltration module
Basic system + UV photooxidation
Basic system + UV photooxidation + ultrafiltration module
Basic system + UV-photooxidation and TOC measurement
Basic system + UV-photooxidation and TOC measurement + ultrafiltration module

GenPure with xCAD wall version:



50131300 standard

50131302 UF

50131317 UV

50131315 UV/UF

50131321 UV-TOC

50131323 UV-TOC/UF

Basic system

Basic system + ultrafiltration module

Basic system + UV-photooxidation

Basic system + UV-photooxidation and ultrafiltration module

Basic system + UV-photooxidation + TOC Measurement

Basic system + UV-photooxidation and TOC Measurement + ultrafiltration module

3 Extent of delivery

Available GenPure with xCAD versions

Intended Use

The GenPure with xCAD 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 GenPure with xCAD 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

Demands on the feedwater

Source	Tap water pretreated by reverse osmosis, ion exchange or distillation.
Silt density index (SDI)	max. 1 for all versions. For water that is not pretreated by reverse osmosis, a prefilter with a 1 µm membrane is recommended.
Feedwater resistance	> 0.5 MΩ _{xc} m
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 to protect the internal filter and, if appropriate, the final filter.
Temperature	2 - 40°C
Pressure	0.1 - 6 bar or 1.45 to 87 PSI

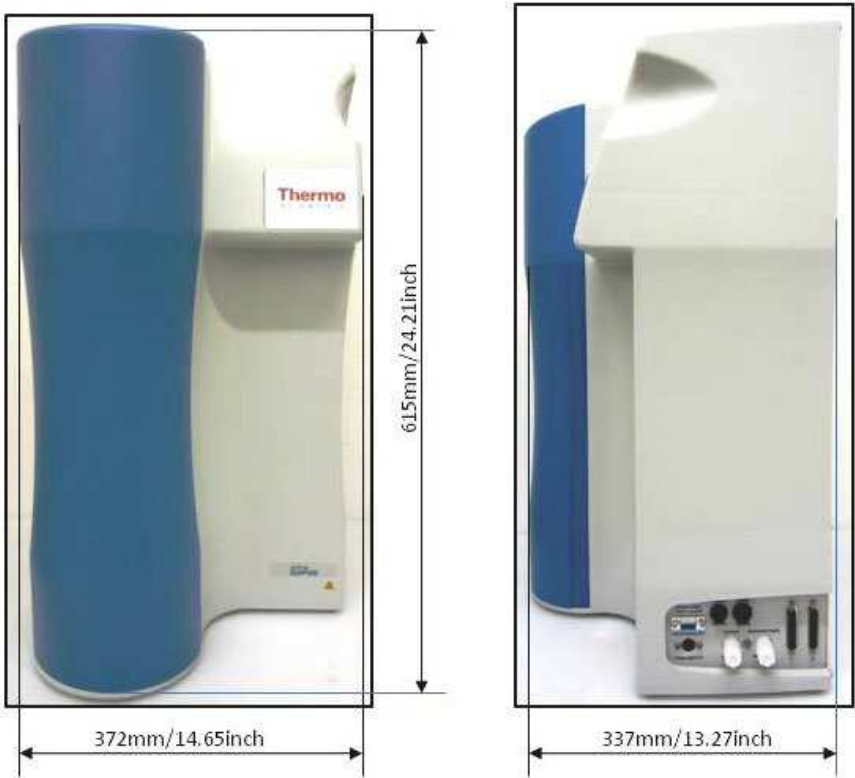
Product water quality

		Standard	UV	UF	UV/UF	UV-TOC	UV-TOC/UF
Resistance (Reference temp. 25 °C)	MΩ_{xc}m at 25 °C	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	< 1	< 1	< 1	< 1	< 1	< 1
Bacterial Endotoxines	EU/ml	--	--	< 0.001*	< 0.001*	--	< 0.001*
Particles	> 0.2 µm	< 1/ml	< 1/ml	< 1/ml	< 1/ml	< 1/ml	< 1/ml

Product water quality							
Flow rate	L/min**	up to 1.5	up to 1.5	up to 1.5	up to 1.5	up to 1.5	up to 1.5
Flow rate with volume control	L/min	1.5	1.5	1.5	1.5	1.5	1.5

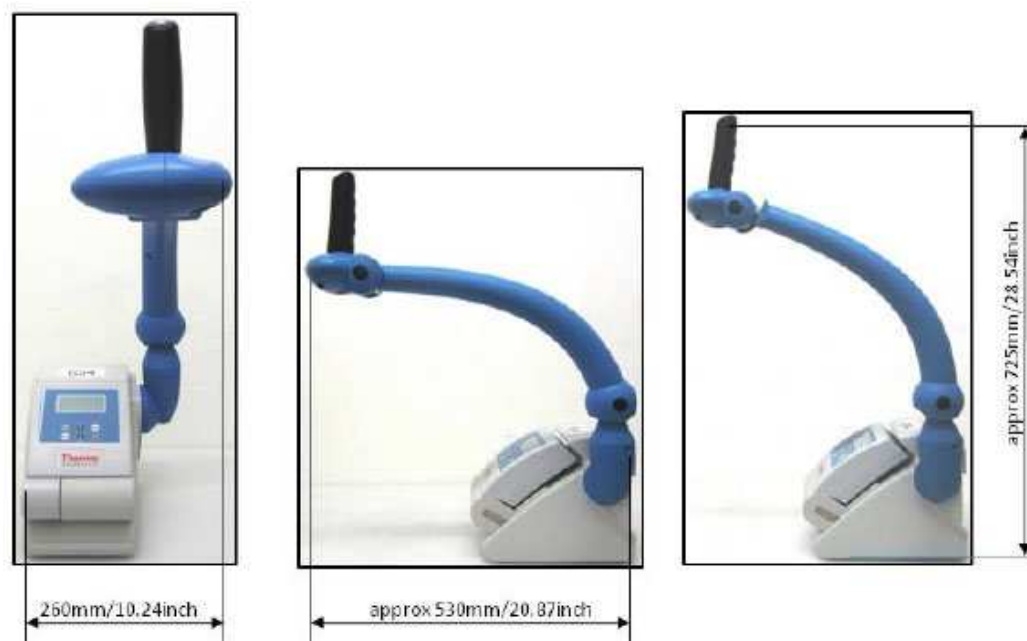
* Depending on feedwater and appropriate disinfection
** Depending on feedwater pressure

Dimension and weight GenPure system		
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)

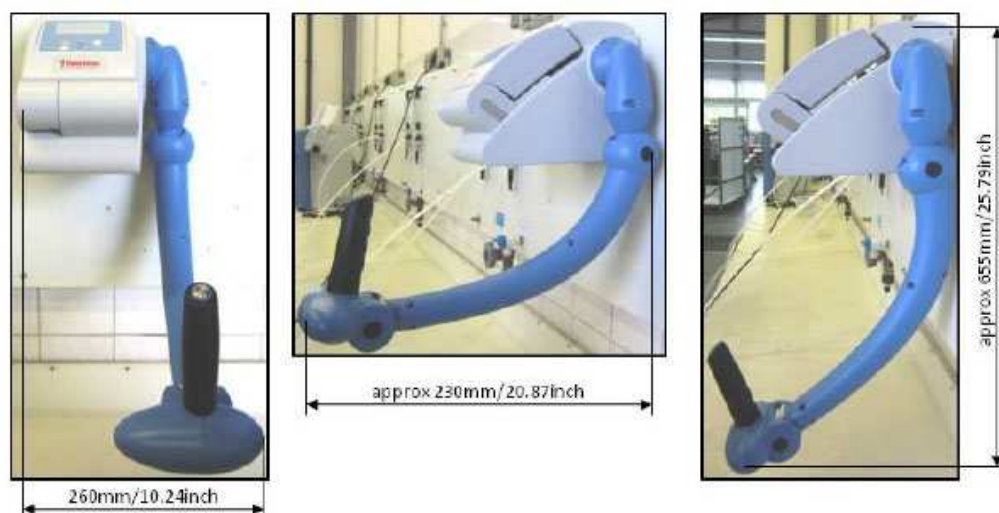


Dimensions and weight xCAD (bench version)

Height	approx. 725 mm	28.54 inch
Width	260 mm	12.24 inch
Depth	approx. 530 mm	20.87 inch
Weight	12 kg	26.46 lbs (dry weight)

**Dimensions and weight xCAD (bench version)**

Height	approx. 655 mm	25.79 inch
Width	260 mm	10.24 inch
Depth	approx. 530 mm	20.87 inch
Weight	5 kg	11.02 lbs (dry weight)



Cell constants, measuring cell

Conductivity, feedwater	0.16 cm ⁻¹
Conductivity, after UV-photooxidation	0.01 cm ⁻¹
Conductivity, ultra pure water	0.01 cm ⁻¹

Connectors for water GenPure

Feedwater	Hose, 8 mm o.d.
Rinse water	Hose, 8 mm o.d.
Ultra pure water/outlet	Hose, 6 mm o.d.
Ultra pure water/recirculation	Hose, 4 mm o.d.

Connectors for water, xCAD

Ultra pure water/inlet	Hose, 6 mm o.d.
Ultra pure water/recirculation	Hose, 4 mm o.d.
Ultra pure water/outlet	R 1/4"
Final filter outlet	Hose, 8 - 10 mm o.d.

Electrical connections / external switched mode power supply

Input voltage	AC 100 – 240 V, 50 – 60 Hz, 5 – 3.8 A
Output voltage	DC 24 V, 3.8 A
System connection	DC 24 V, 80 W
Serial interface	RS 232
Protection Class	Class II (external SMPS certified as Class I)

Electrical connections, xCAD

2x SUB-D socket	25 pin
-----------------	--------

Airborne sound emission

Sound-pressure level	49 db(A)
----------------------	----------

Ambient conditions		
Usage		Indoor rooms
Height		Up to 2000 m
Temperature range		From 5 °C to 40 °C
Relative humidity		Maximum relative humidity 80 % at temperatures of up to 31 °C, linearly decreasing to 50 % relative humidity at 40 °C
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).
<div style="text-align: center;"> <div style="background-color: #0056b3; color: white; padding: 5px; display: inline-block;">NOTE</div> <p>The rated level of transient overvoltage is the withstand impulse voltage acc. to overvoltage category II of IEC 60364-4-443</p> </div>		
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 fibre
UV lamp		High purity quartz
UV housing		Stainless steel
Ultrapure cartridge		PP = polyethylene
UF housing		PC = polycarbonate
Rinsing solenoid valve		PA = polyamid
Dispensing valve		PVDF = polivinyldenfluorid
Conductivity measuring cell		POM = polyoxymethylen, stainless steel
Distributor block		POM = polyoxymethylen
Connectors		POM = polyoxymethylen
Hoses		PE = polyethylene
O-rings		EPDM = ethylen-propylen-diene-rubber

The installation area

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).



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.
- Level standing surface.
- 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 specification](#)“ on [page 18](#)).
- A floor drain with a nominal out diameter of 63mm or 2.48 inch (DN 50 pipe) drain pipe.
- Free run off to drain.
When no floor drain is available, install a water watcher to protect against water damage (only available for EU).



Unrestricted gravity flow to drain must be ensured!

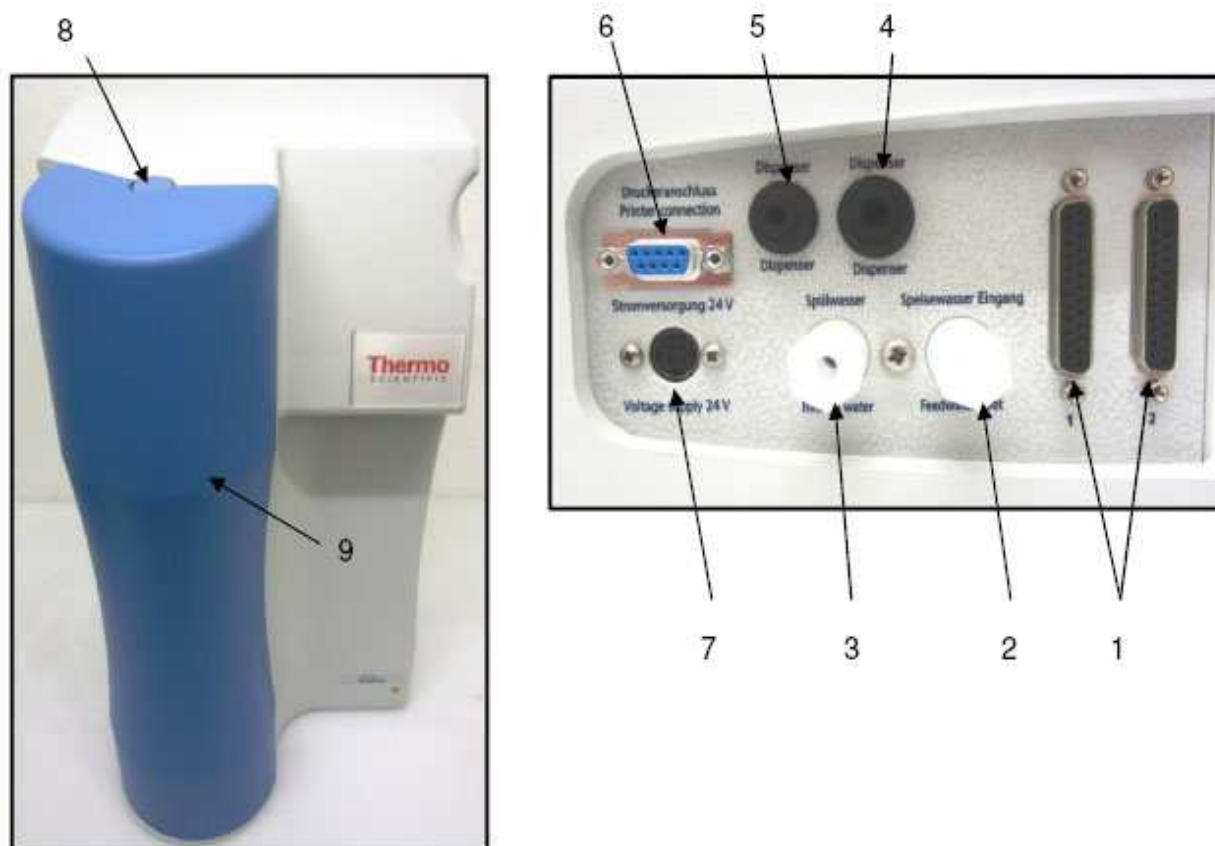
- An electrical socket appropriate for the system (see „[Technical specification](#)“ on [page 20](#)).
- Sufficient working room all around the system (approx 30 cm / 11.81 inch, for replacing filters etc.).
- Easy access for operation and control of the system.
- Water pre treated such as DI, RO or distillation water connection with 3/4 NPT male thread and customer supplied shutoff valve.

Installation

Contents

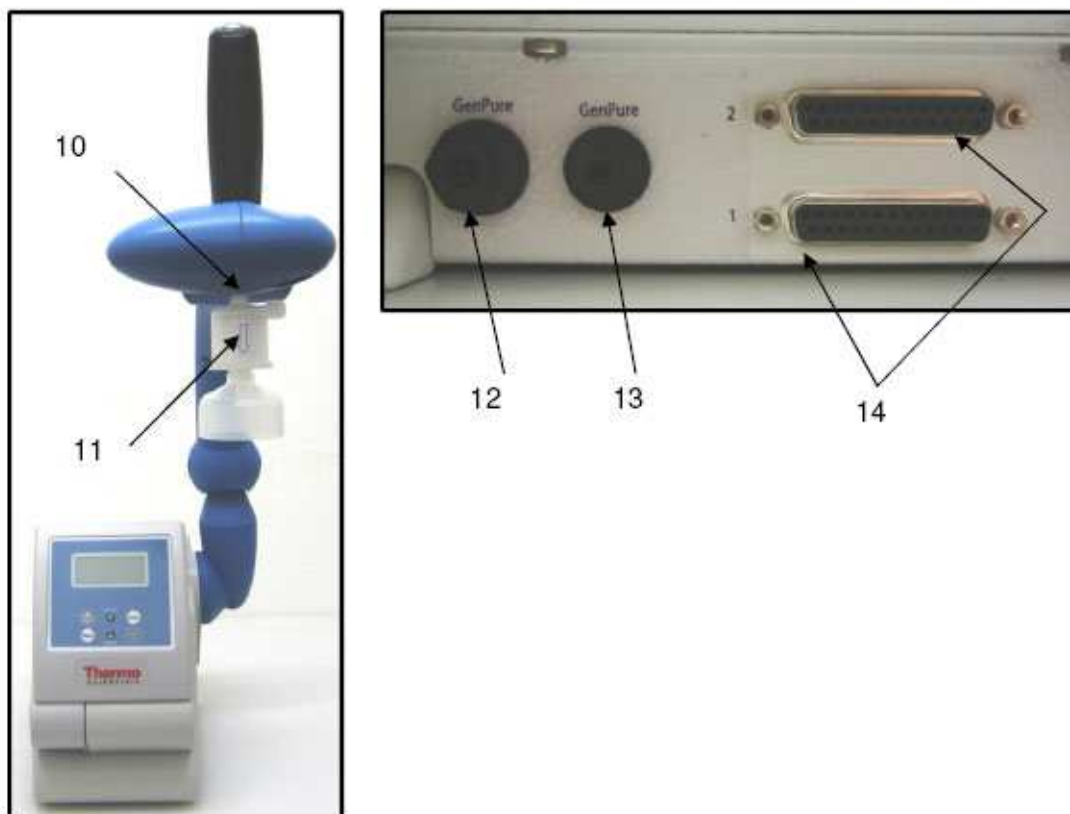
- “Installation of GenPure with xCAD system, bench version” on page 28
- “Installation of GenPure with xCAD system, wall version” on page 32
- “Wall mounting GenPure system” on page 37
- “Mounting the power pack (voltage supply)” on page 39
- “Installation examples” on page 41

Connectors GenPure system



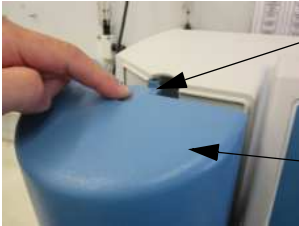

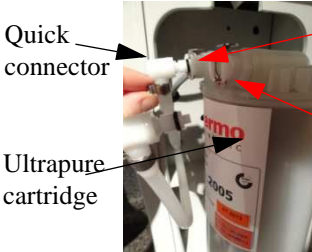
1. Connectors for 25 pin sockets to xCAD (system control)
2. Feedwater connector, 8 mm o.d or 0.31"
3. Rinse water connector, 8 mm o.d 0.31"
4. Ultrapure water connector, 6 mm o.d or 0.24" (to xCAD)
5. Ultrapure water connector, 4 mm o.d 0.16" (to xCAD)
6. Optional printer connection
7. Power supply connector 24 V DC
8. Push button for releasing the cartridge cover
9. Cartridge cover


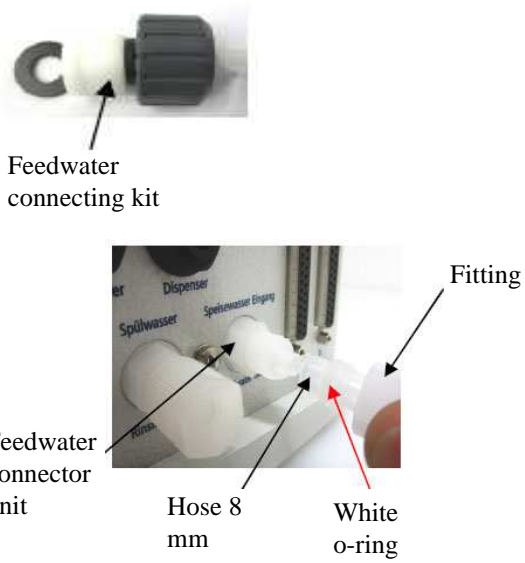
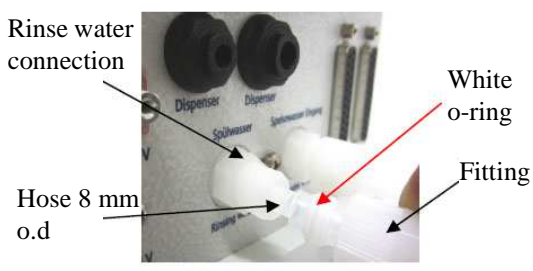
Connectors xCAD

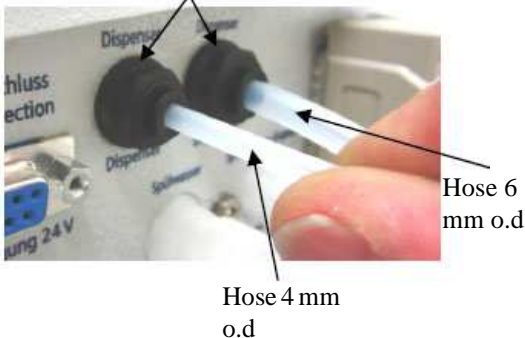
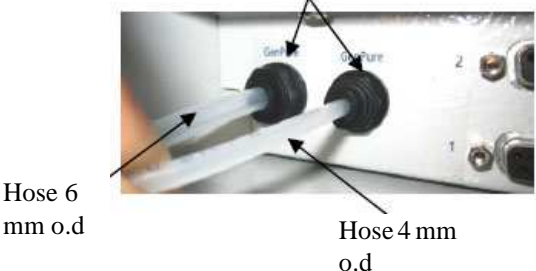
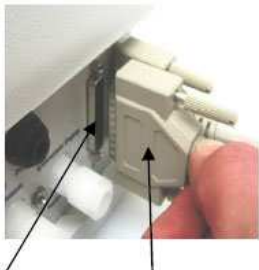

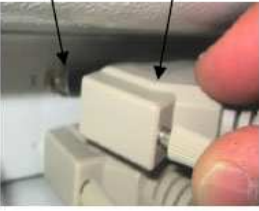






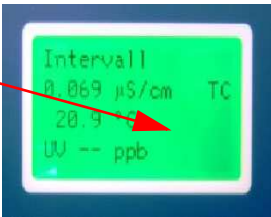
- 10. Dispensing valve outlet, R 1/4" female thread
- 11. Final filter 0.2 μm
- 12. Ultrapure water connector, 6 mm o.d or 0.24" (to GenPure)
- 13. Ultrapure water connector, 4 mm o.d or 0.16" (to GenPure)
- 14. Connector for 25 pin sockets to GenPure (system control)

Installation of GenPure with xCAD system, bench version

Step	Action	Figure
1	Either place the GenPure system with the xCAD bench version on the intended surface or hang it on a wall. For wall mounting the GenPure system using the included wall mounting hardware.	<div data-bbox="1173 376 1460 450" data-label="Text"> <p>NOTE</p> </div> <p>See under chapter “Wall mounting GenPure system” on page 37.</p>
2	Release the cartridge cover by pressing the push button.	 <p>Push-button</p> <p>Cartridge cover</p>
3	Remove the two stoppers from the new ultrapure cartridge and fit the cartridge into the system.	 <p>Stoppers</p>
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 connector</p> <p>Outlet</p> <p>Inlet</p> <p>Ultrapure cartridge</p>

Step	Action	Figure
5	<p>.Mount the feedwater connecting kit together and connect it to the feedwater inlet line. Connect the other end of the hose to the feedwater connector of the system by unscrewing the fitting. After this put the hose through the fitting and mount the white O-ring on it. Screw the fitting back to the system.</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> <p>Fitting</p> <p>Feedwater connector unit</p> <p>Hose 8 mm</p> <p>White o-ring</p>
6	<p>Connect the 8mm o.d hose to the rinse water connection of the system (see step 5) and make a gravity fall (pressureless) connection from the system to the floor drain. The drain to the sewer must be max. 1m (1.09 yards) above the rinsing water connector of the unit.</p>	 <p>Rinse water connection</p> <p>Hose 8 mm o.d</p> <p>White o-ring</p> <p>Fitting</p>

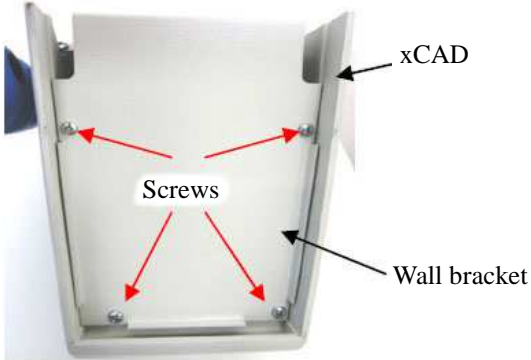
Step	Action	Figure
7	<p>NOTE</p> <p>Before you begin with step a guide the two hoses 6 and 4 mm o.d or 0.16 and 0.24 inch through the cable spiral.</p> <ol style="list-style-type: none"> Connect the two ends of the hoses 6 and 4 mm o.d or 0.16 and 0.24 inch of the cable spiral into the ultrapure water connectors of the GenPure system. The other ends of the hoses 6 and 4 mm o.d or 0.16 and 0.24 inch you should connect to the connectors of the xCAD. 	<p>a) Connectors GenPure</p>  <p>b) Connectors xCAD</p> 
8	<ol style="list-style-type: none"> Plug each of the cables with 25 pin socket in the sockets of the GenPure system and screw them tight. Plug in the two 25 pin system control cable onto the xCAD and fix them with the screws <p>NOTE</p> <p>In order to recognise the correct connectors of the cables on the xCAD of system and the xCAD they are marked with numbers 1 and 2 (see red arrows).</p>	<p>a)</p>  <p>25 pin connectors</p>  <p>b)</p>  <p>25 pin connectors</p> 

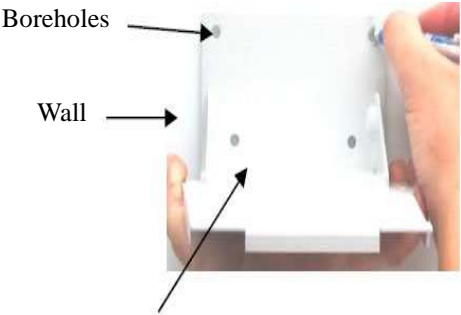
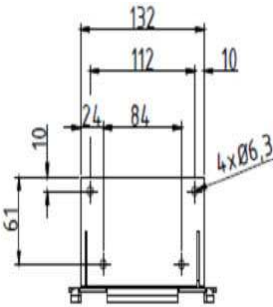
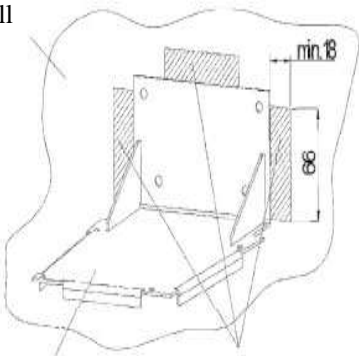
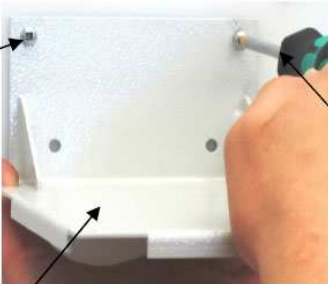
Step	Action	Figure
9	Screw the final filter in counter clockwise direction (see red arrow in the picture) into the 1/4" female thread of the xCAD dispensing valve.	 <p>1/4" female thread connection</p> <p>Final filter 0.2 μm</p>
10	Assemble the power pack and make the voltage connection to the GenPure system.	<div data-bbox="906 683 1193 759" data-label="Text">NOTE</div> <p>See under chapter “Mounting the power pack (voltage supply)” on page 39.</p>
11	<p>If applicable use the RS232 connector (6) to connect the optional data printer.</p> <p>Open the feedwater supply tap.</p> <div data-bbox="325 1030 604 1104" data-label="Image"> <p>CAUTION</p> </div> <p>Only feedwater that has been pretreated by reverse osmosis, ion exchange or distillation is to be used.</p>	   <p>Feedwater supply</p>

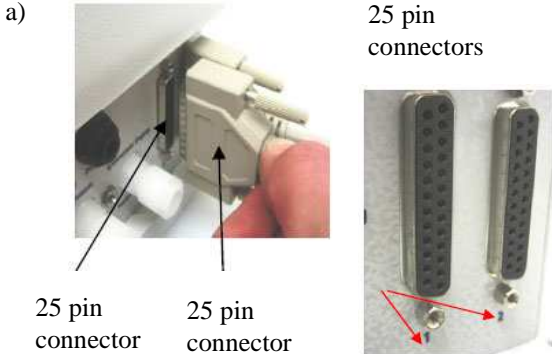
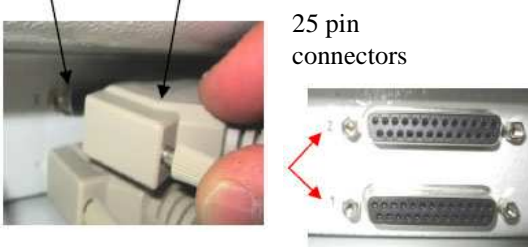
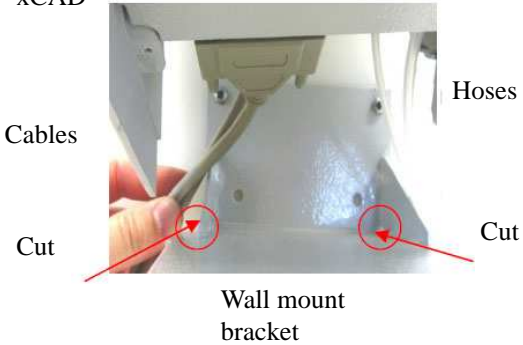
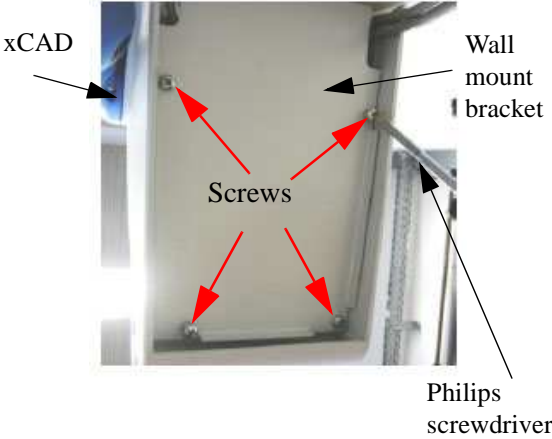
Installation of GenPure with xCAD system, wall version

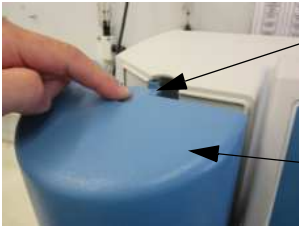

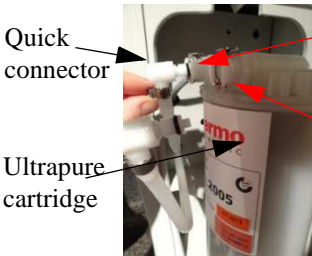

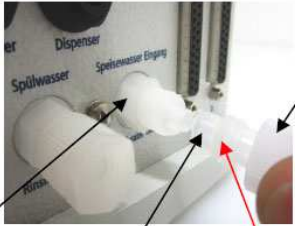



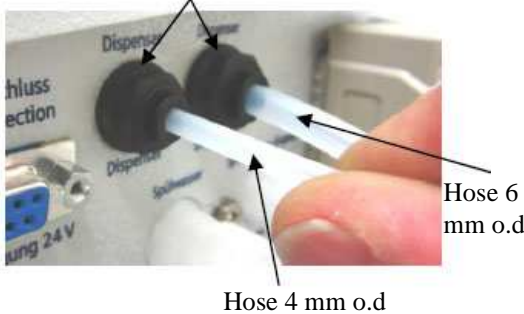
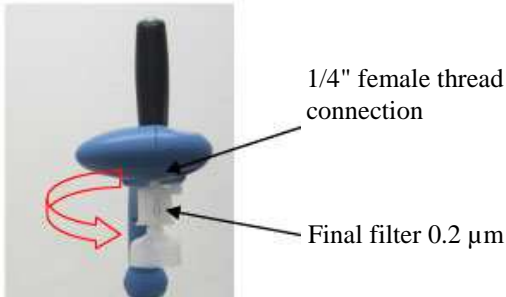


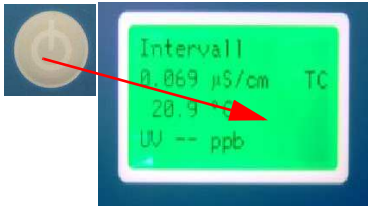
Before hanging the xCAD onto a wall make sure that the wall can support the weight of the system once it's full of water.

Step	Action	Figure
1	Either place the GenPure system on the intended surface or hang it on a wall. For wall mounting the GenPure system using the included wall mounting hardware.	<div data-bbox="906 607 1193 683" data-label="Section-Header"> <h3>NOTE</h3> </div> <p>Lift and carried out the xCAD wall version by two people. It is easier to work and mount it onto a wall.</p>
2	To wall mount the xCAD wall version unscrew the 4 screws (see red arrows in the picture) of the underside from the xCAD and remove the wall mount bracket.	

Step	Action	Figure
3	<p>a. Hold the wall mount bracket at the desired position on the wall and mark the four boreholes for fixing the wall mount bracket.</p> <p>Then use a 6 mm or 1/4" twist drill to make the holes and put in the four S6 dowels which are supplied in the assembly kit.</p> <p>NOTE</p> <p>If you are want to take the hoses and cables out of back the wall look at the pictures 1) and 2) and step b). When it is not wish going to step 4.</p> <p>b. Refer to dimensions on picture 1) and 2) to make the necessary wall cuts needed if you want to push the 0.31" (8mm) o.d hoses and cable out through the wall behind the xCAD.</p>	<p>a)</p>  <p>1)</p> <p>Dimensions boreholes of wall mount bracket</p>  <p>2)</p> <p>Wall</p>  <p>Wall mount</p> <p>Possible wall cut- outs for cable and hose taken out at back</p>
4	<p>Attach the wall mount bracket to the wall by screwing in the 4 supplied screws with a philips screw driver into the wall where you have put in the plugs before.</p>	 <p>Screw</p> <p>Philips screw-driver</p> <p>Wall mount bracket</p>

Step	Action	Figure
5	<p>NOTE</p> <p>Before you begin with step 5 guide the two hoses 6 and 4 mm o.d or 0.16 and 0.24 inch through the cable spiral.</p> <ol style="list-style-type: none"> C. Plug each of the cables with 25 pin socket in the sockets of the GenPure system and fix them with the screws. Plug in the two 25 pin system control cable onto the system and fix them with the screws. <p>NOTE</p> <p>In order to recognise the correct connectors of the cables on the xCAD of system and the xCAD are they marked with numbers 1 and 2 (see red arrows).</p>	<p>a)</p>  <p>25 pin connectors</p> <p>25 pin connector 25 pin connector</p> <p>b)</p>  <p>25 pin connectors</p>
6	<p>Place the xCAD wall version onto the mounted wall mount bracket. There are two cuts on the bracket (see red arrows) the cables and hoses for guidance.</p> <p>NOTE</p> <p>When you have made the possible wall cuts (see step 3) plug the cables and hoses throughout the wall.</p>	<p>xCAD</p>  <p>Cables</p> <p>Hoses</p> <p>Cut Cut</p> <p>Wall mount bracket</p>
7	<p>Screw in the 4 screws (see red arrows) which you unscrewed in step 2 to attach the xCAD Client on the wall mount bracket.</p>	 <p>xCAD</p> <p>Wall mount bracket</p> <p>Screws</p> <p>Philips screwdriver</p>

Step	Action	Figure
8	Release the cartridge cover by pressing the push button.	 <p>Push-button</p> <p>Cartridge cover</p>
9	Remove the two stoppers from the new ultrapure cartridge and insert the cartridge into the system.	 <p>Stoppers</p>
10	<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 connector</p> <p>Outlet</p> <p>Inlet</p> <p>Ultrapure cartridge</p>
11	<p>Mount the feedwater connecting kit together and connect it to the feedwater sample. Connect the other end of the hose from the feedwater connecting kit to the feedwater connector of the system by unscrewing the fitting. After this stick the hose through the fitting and mount the white o-ring on it. Screw the fitting back to the system</p>	 <p>Feedwater connecting kit</p>
	<p>CAUTION</p> <p>Only feedwater that has been pretreated by reverse osmosis, ion exchange or distillation is to be used.</p>	 <p>Fitting</p> <p>Feedwater connector unit</p> <p>Hose 8 mm</p> <p>White o-ring</p>

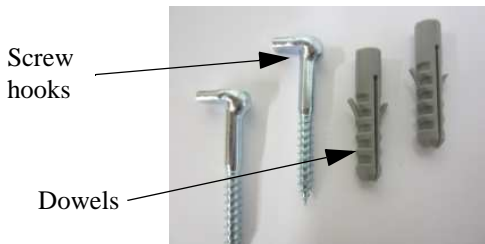

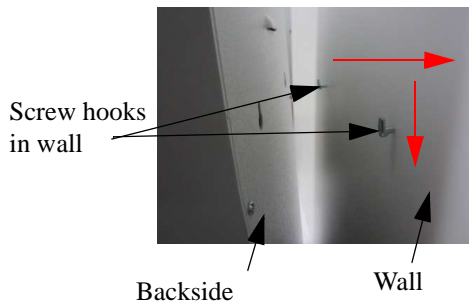
Step	Action	Figure
12	Connect the 8mm o.d (0.31")hose to the rinse water connection of the system (see step 5) and make a gravity fall (pressureless) connection from the system to the floor drain. The drain to the sewer must be max. 1m(1.07 yards) above the rinsing water connector of the unit.	
13	Connect the two other ends of the hoses 6 and 4 mm o.d or 0.16 and 0.24 inch of the cable spiral which you have mounted in step 5 to the xCAD into the ultrapure water connectors of the GenPure system.	
15	Screw the final filter in counter clockwise direction (see red arrow in the picture) into the 1/4" female thread of the xCAD dispensing valve.	
16	Assemble the power pack and make the voltage connection to the GenPure system.	<div data-bbox="906 1435 1193 1509" data-label="Text">NOTE</div> <p>See under chapter “Mounting the power pack (voltage supply)” on page 39.</p>
17	<p>If applicable use the RS232 connector (6) to connect the optional data printer.</p> <p>Open the feedwater supply tap.</p> <div data-bbox="328 1776 593 1845" data-label="Image">  </div> <p>Only feedwater that has been pretreated by reverse osmosis, ion exchange or distillation is to be used.</p>	 

Wall mounting GenPure system

NOTE

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. Then use a twist drill (8mm 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 that are supplied in the assembly kit in the holes. Screw the 5.2 x 50mm screw hooks that are also supplied in the assembly kit into the dowels.	
3	Lift the GenPure System and hang the back side of it onto the screw hooks.  Lifting and carrying the GenPure system should be completed by 2 people.	

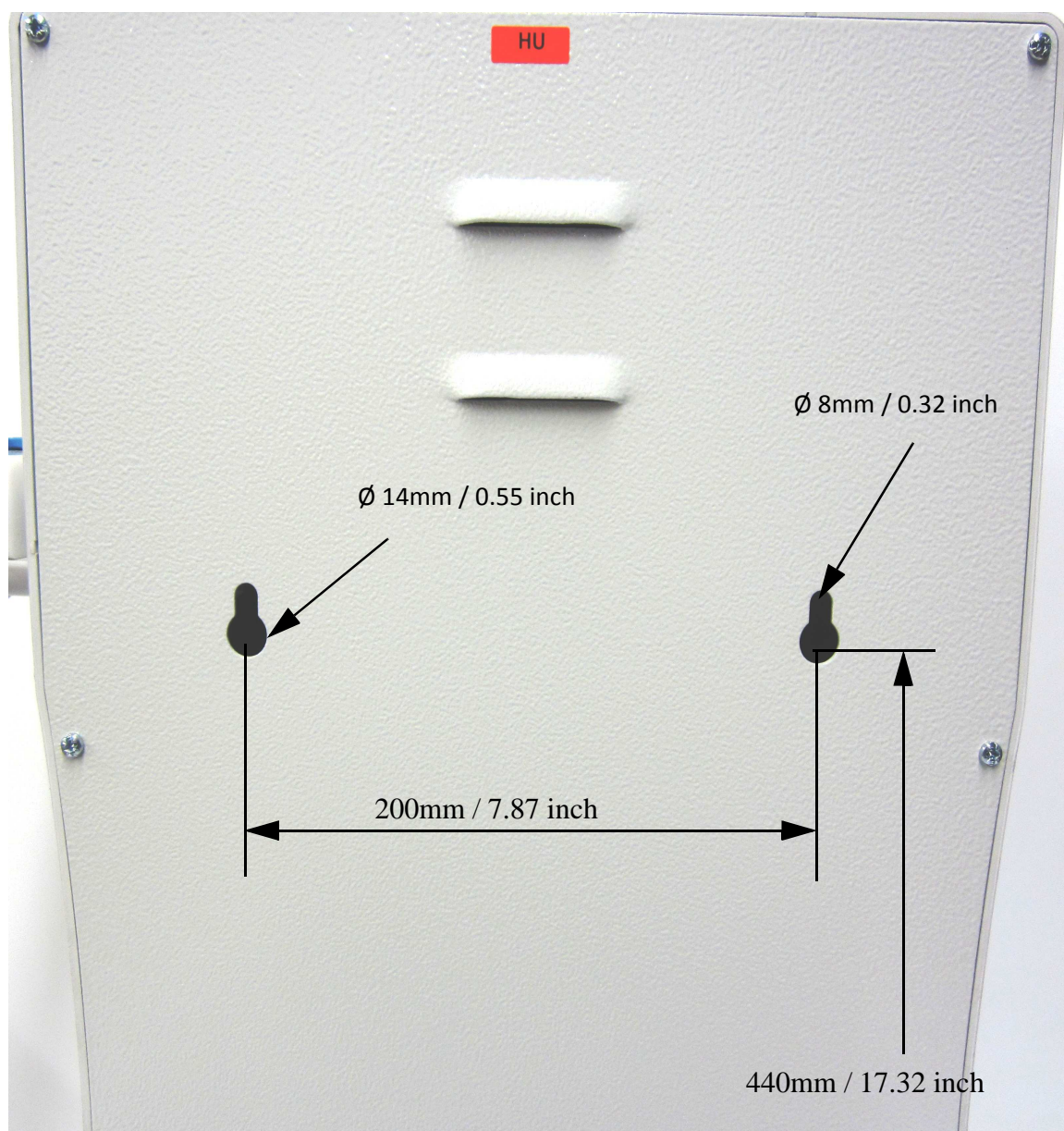
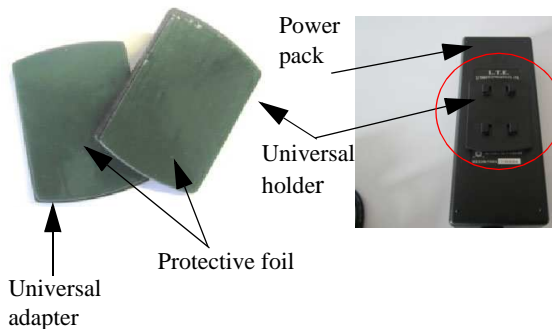
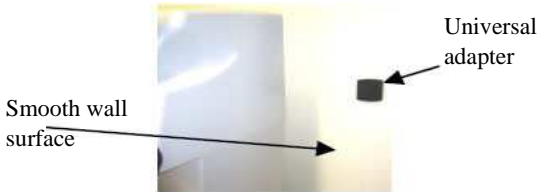
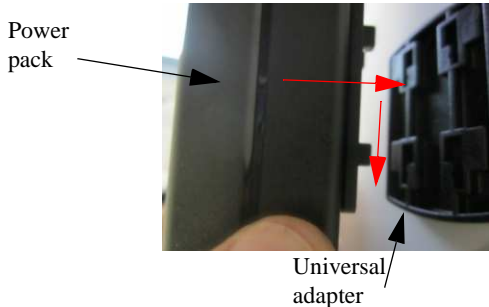
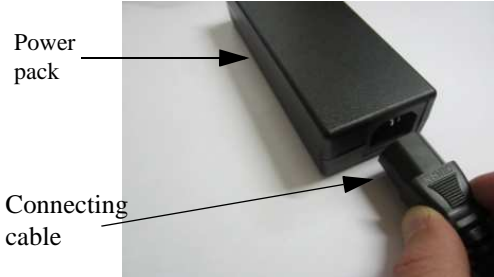



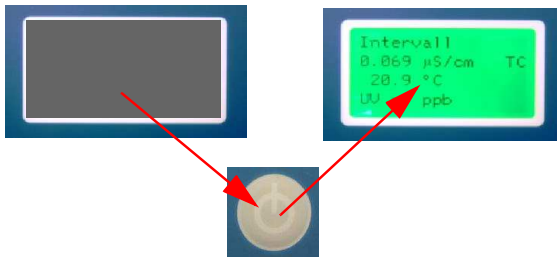
Figure 1. Holes for wall mounting

Mounting the power pack (voltage supply)

NOTE

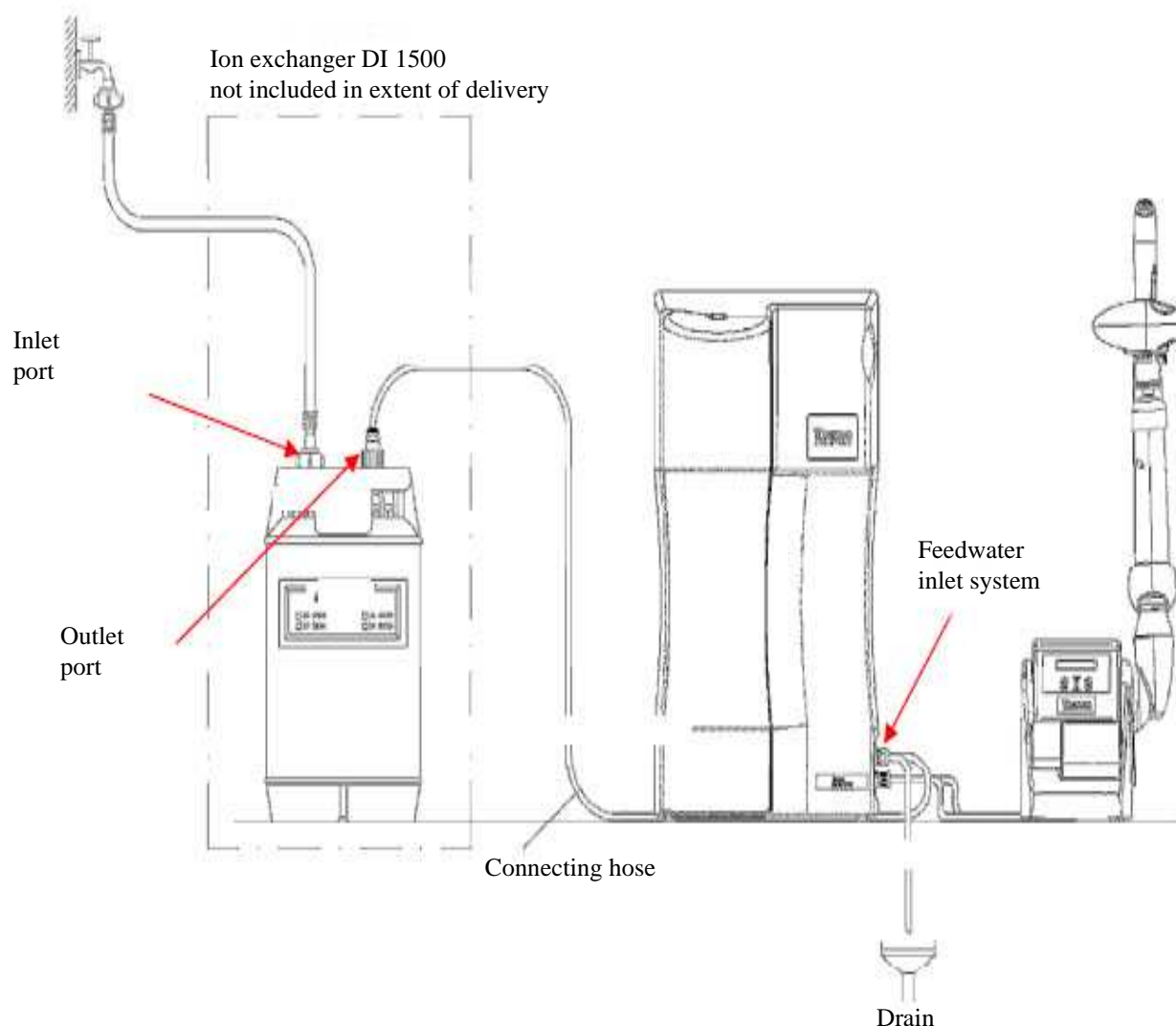
Whenever possible, mount the power pack on the wall to the left or right of the ultra pure water system where it is freely accessible and will not come in contact with water for get wet.

Step	Action	Figure
1	<p>NOTE</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 above figure next to this text.</p>	 <p>Power pack</p> <p>Universal holder</p> <p>Universal adapter</p> <p>Protective foil</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>	 <p>Universal adapter</p> <p>Smooth wall surface</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>NOTE</p> <p>The removable line cord must be shown to the bottom.</p>	 <p>Power pack</p> <p>Universal adapter</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>	 <p>Power pack</p> <p>Connecting cable</p>

Step	Action	Figure
5	Connect the power pack to the ultrapure water system (24V 4-pin power supply connector, connector 8) and to an earthed 100 - 250V, 50/60Hz socket.	 <p>Power supply connector</p>
6	Put the system on. The system is now ready for use.	

Installation examples

Connection to an 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.

Flow charts

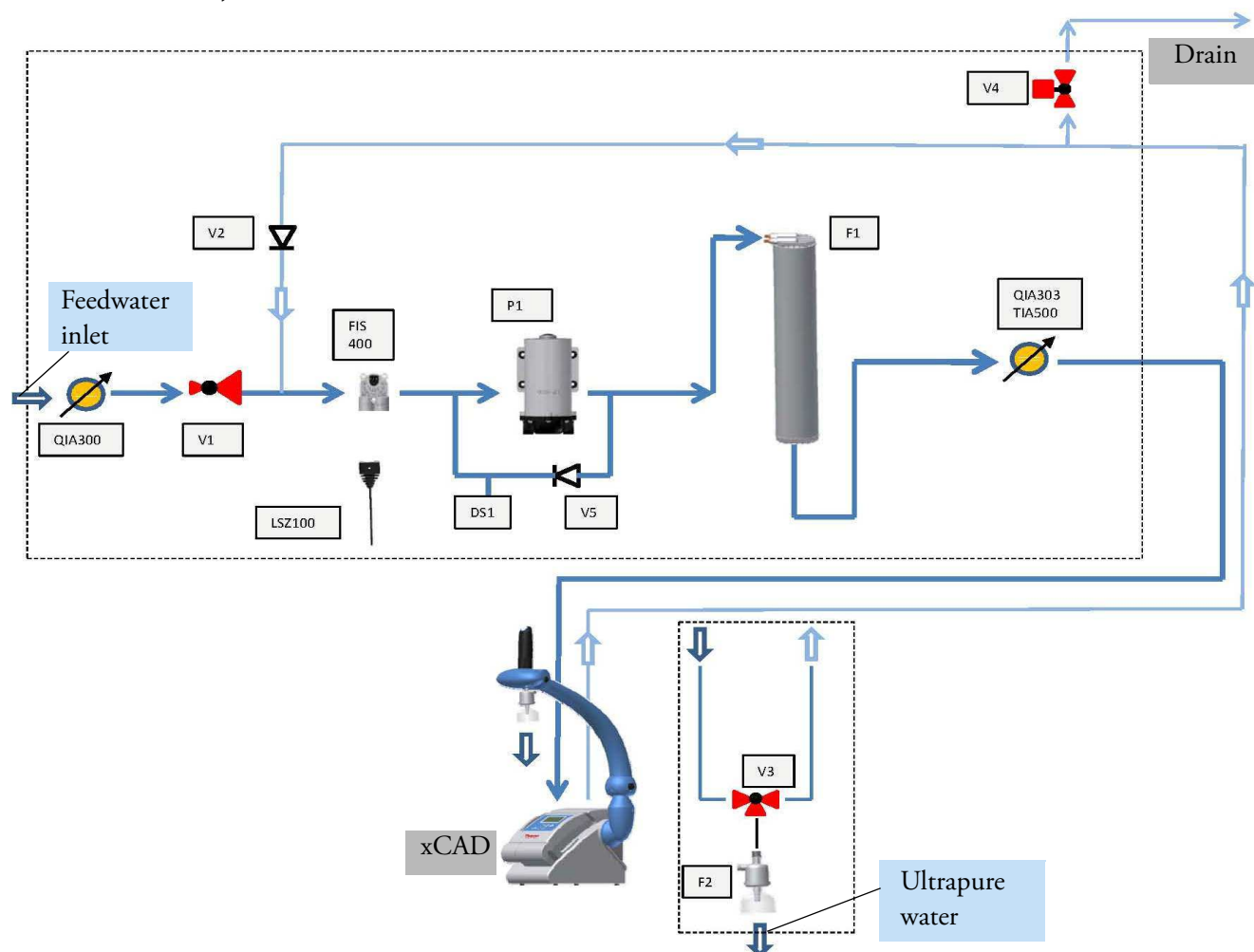
Contents

- “Flow chart, GenPure with xCAD standard” on page 43
- “Flow chart, GenPure UV with xCAD” on page 44
- “Flow chart for GenPure UF” on page 45
- “Flow chart, GenPure UV/UF with xCAD” on page 46
- “Flow chart, GenPure UV-TOC with xCAD” on page 47
- “Flow chart, GenPure UV-TOC/UF with xCAD” on page 48

8 Flow charts

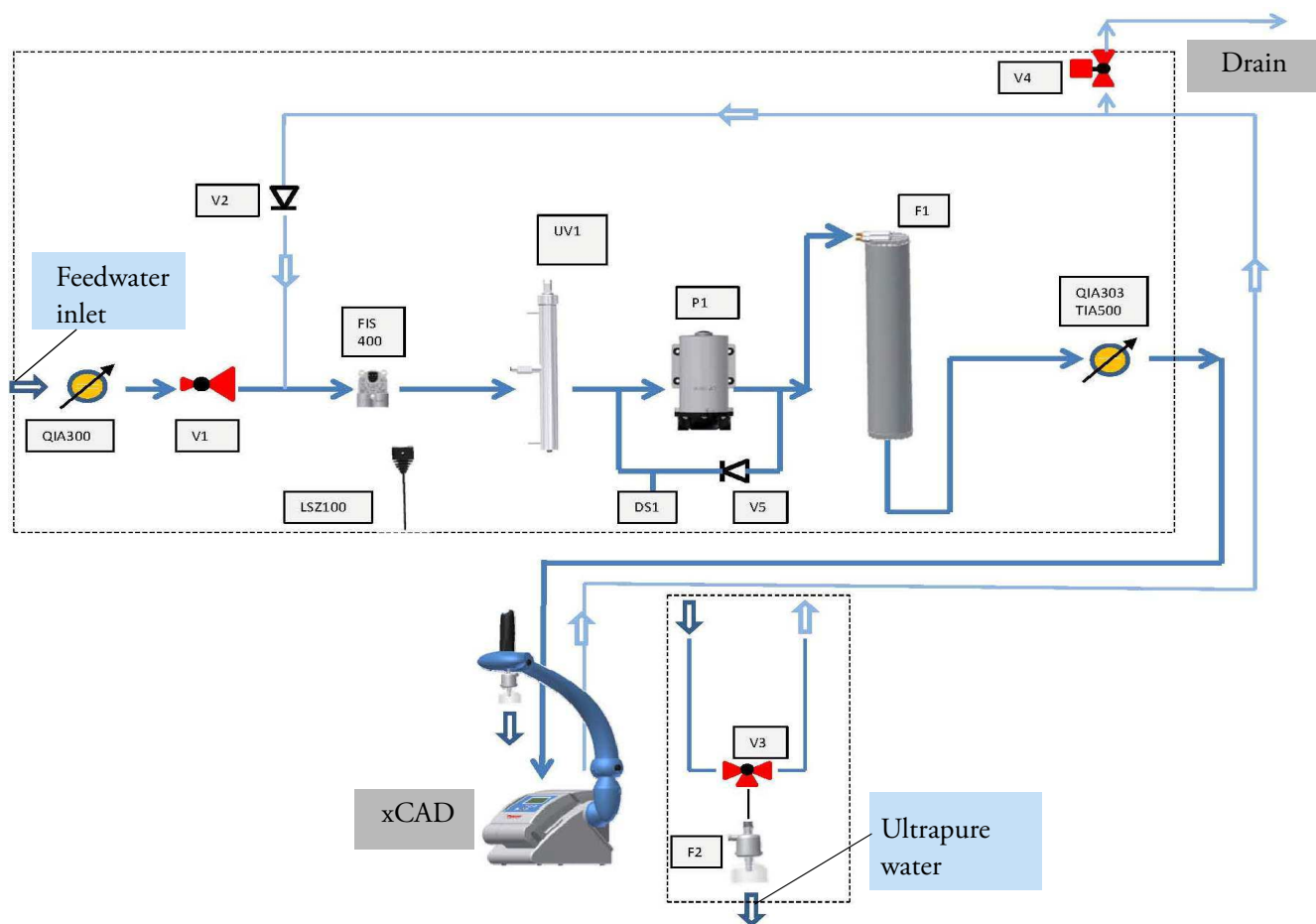
Flow chart, GenPure with xCAD standard

Flow chart, GenPure with xCAD standard



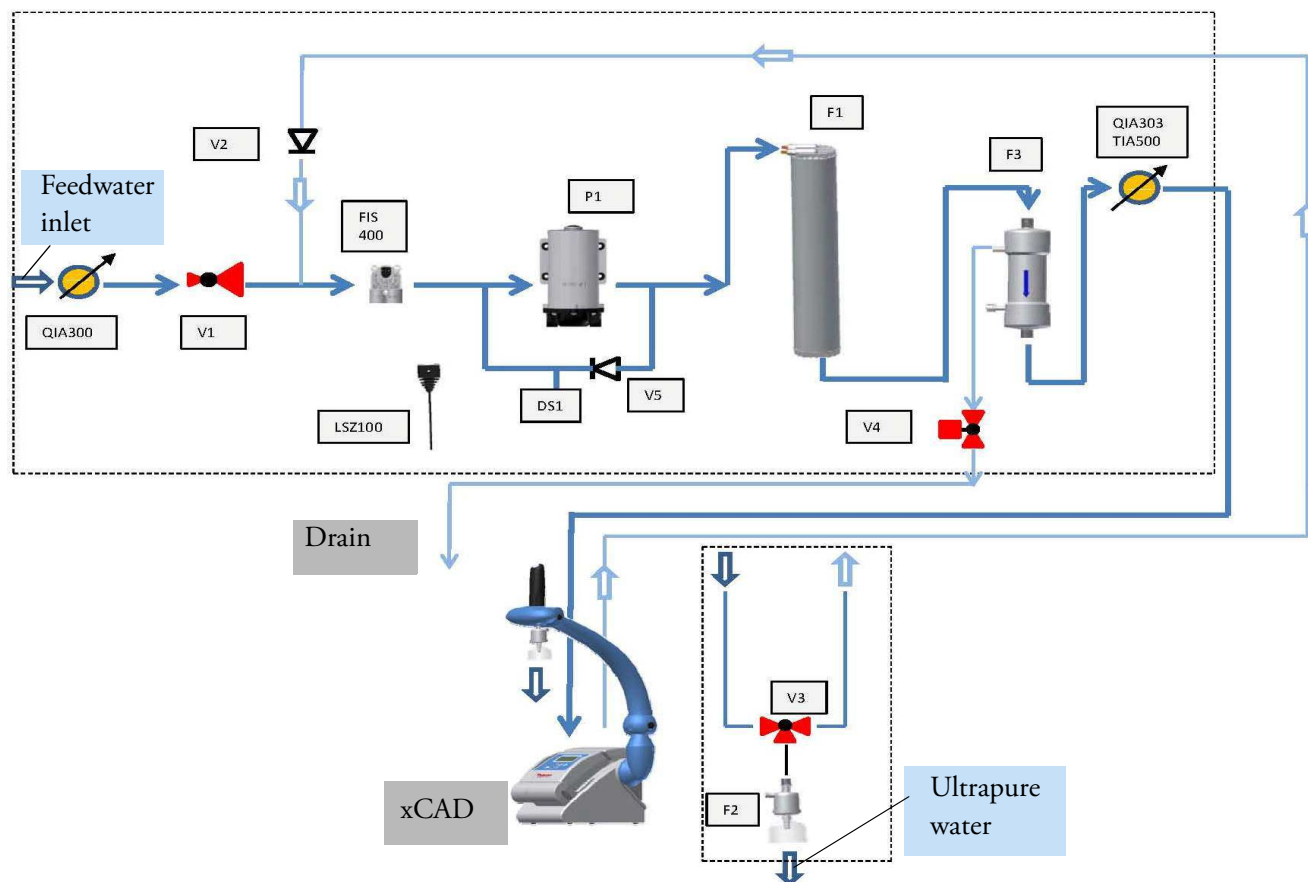
- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- P1 Circulation pump
- FIS400 Digital flowmeter
- QIA 300 Conductivity, feedwater
- QIA 301 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

Flow chart, GenPure UV with xCAD



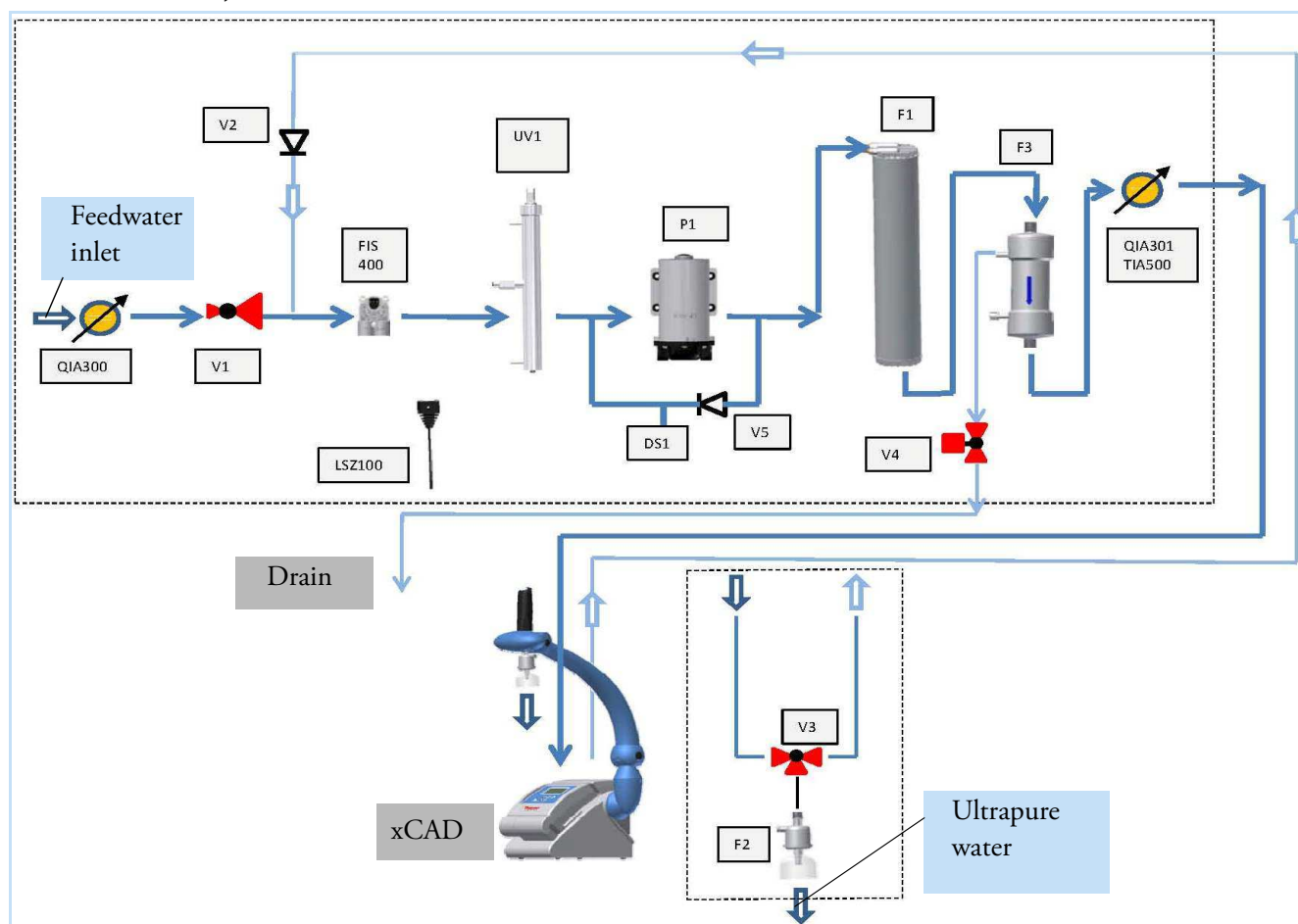
- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- P1 Circulation pump
- UV1 UV-photooxidation
- FIS400 Digital flowmeter
- QIA 300 Conductivity feedwater
- QIA 301 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

Flow chart for GenPure UF



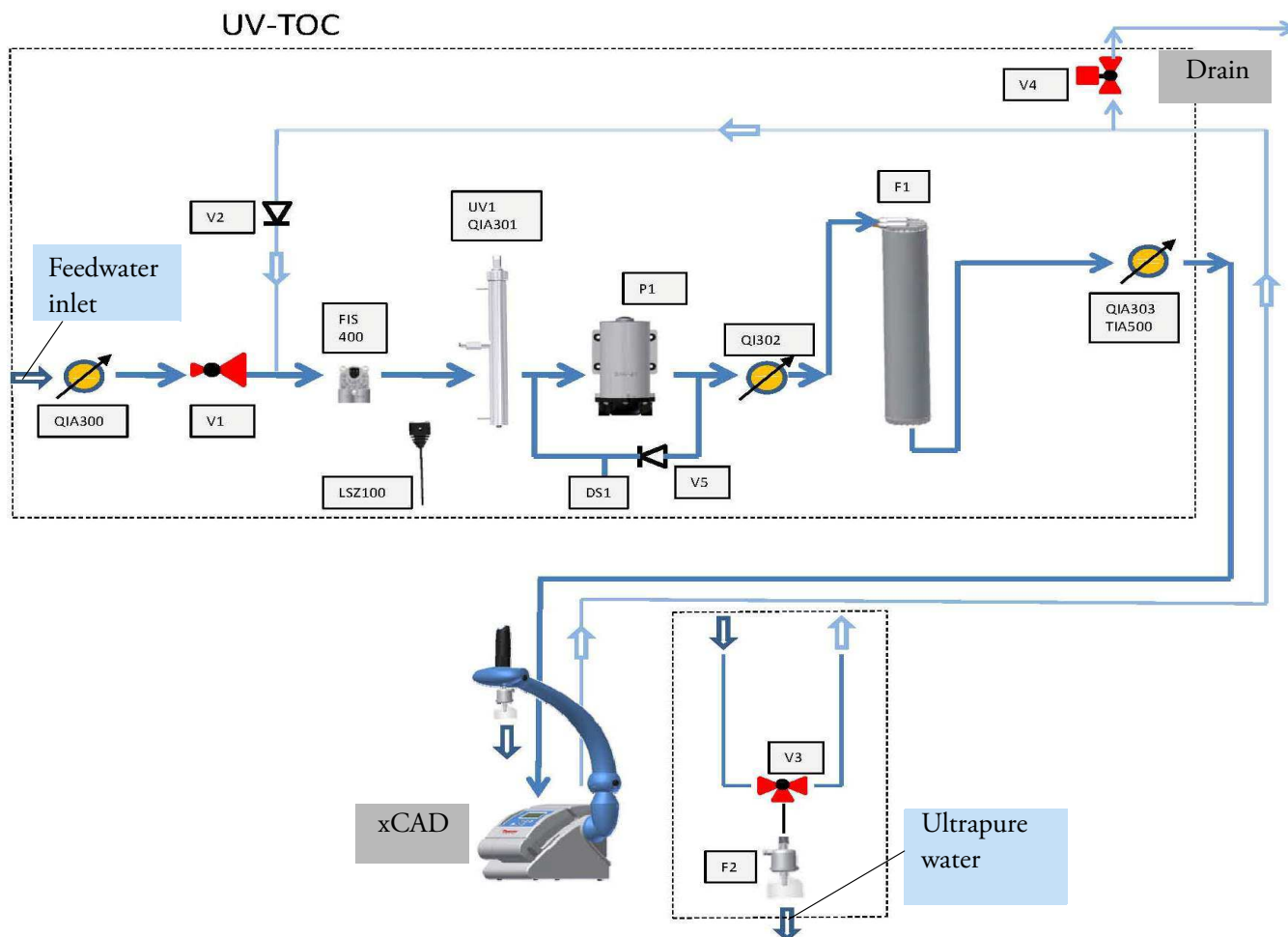
- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- F3 Ultrafiltration module
- P1 Circulation pump
- FIS400 Digital flowmeter
- QIA 300 Conductivity, feedwater
- QIA 301 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

Flow chart, GenPure UV/UF with xCAD



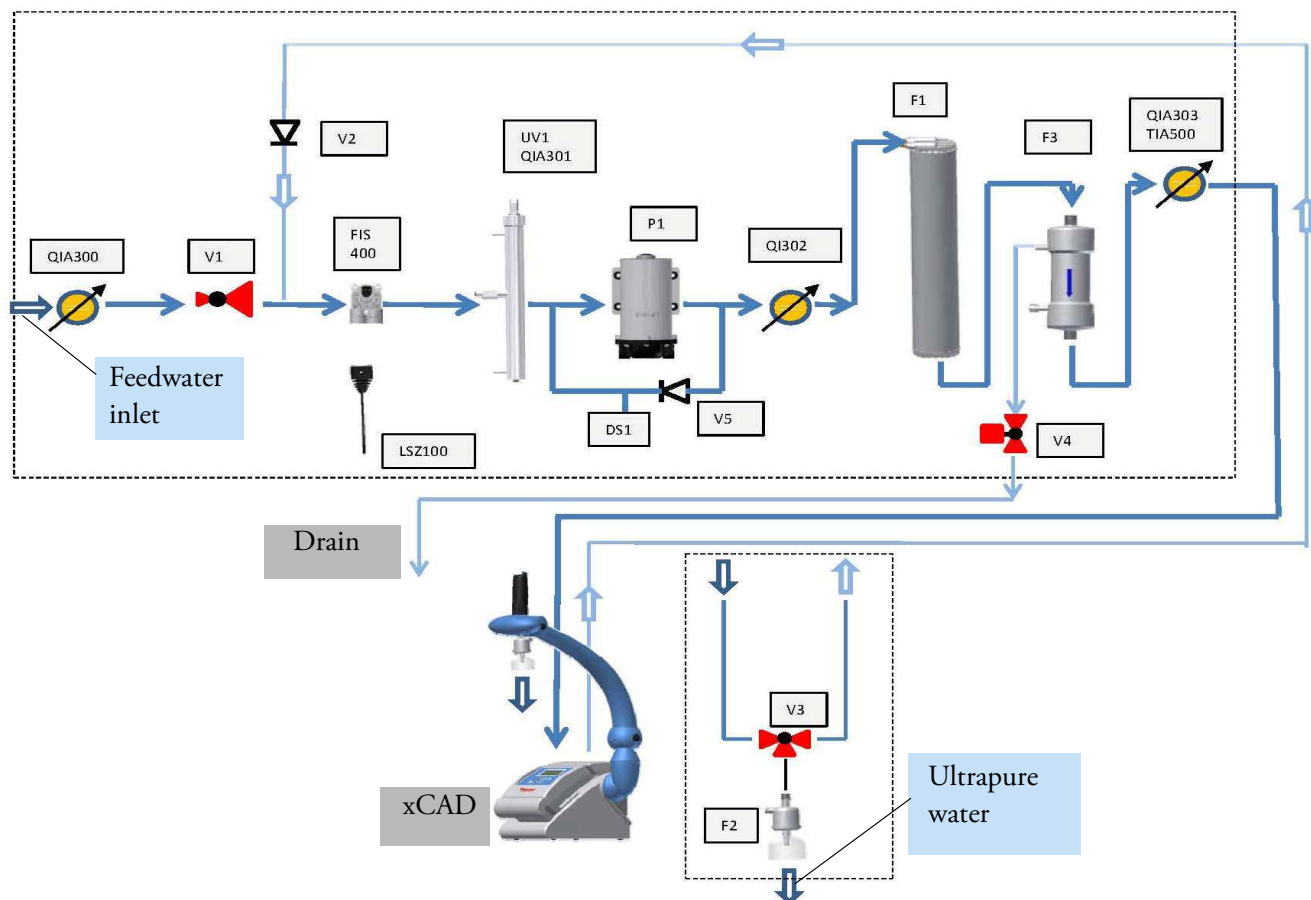
- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- F3 Ultrafiltration module
- P1 Circulation pump
- UV1 UV-photooxidation
- FIS 400 Digital flowmeter
- QIA 300 Conductivity, feedwater
- QIA 301 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

Flow chart, GenPure UV-TOC with xCAD



- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- P1 Circulation pump
- UV1 UV-photoreactor
- FIS400 Digital flowmeter
- QIA 300 Conductivity, feedwater
- QIA 301 UV-Intensity
- QI 302 Conductivity, TOC measurement
- QIA 303 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

Flow chart, GenPure UV-TOC/UF with xCAD



- DS1 Dosing orifice
- F1 Ultrapure cartridge
- F2 Final filter
- F3 Ultrafiltration module
- P1 Circulation pump
- UV1 UV-photooxidation
- FIS400 Digital flowmeter
- QIA 300 Conductivity, feedwater
- QIA 301 UV-Intensity
- QI 302 Conductivity, TOC measurement
- QIA 303 Conductivity, ultra pure water
- TIA 500 Temperature sensor
- V1 Pressure reducer
- V2 Check valve
- V3 Solenoid valve
- V4 Rinsing solenoid valve
- V5 Check valve 1 bar

8 Flow charts

Flow chart, GenPure UV-TOC/UF with xCAD

How the system function

NOTE

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 a end 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 a 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.

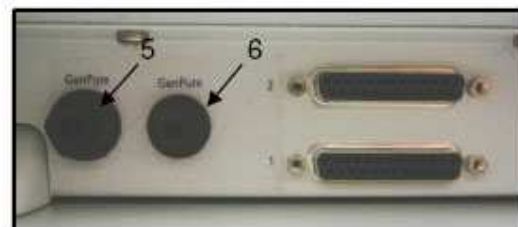
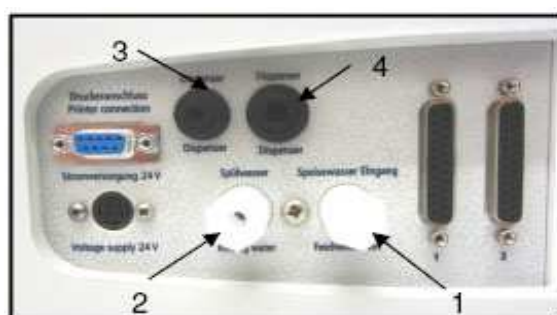
Putting system into operation

NOTE

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 connection system 8 mm o.d or 0.31"
2. Rinse water connection system 8 mm o.d or 0.31"
3. Ultrapure water connection system 4 mm o.d or 0.16"
4. Ultrapure water connection system 6 mm o.d or 0.24"
5. Ultrapure water connection xCAD 6 mm o.d or 0.24"
6. Ultrapure water connection xCAD 4 mm o.d or 0.16"



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

NOTE

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.



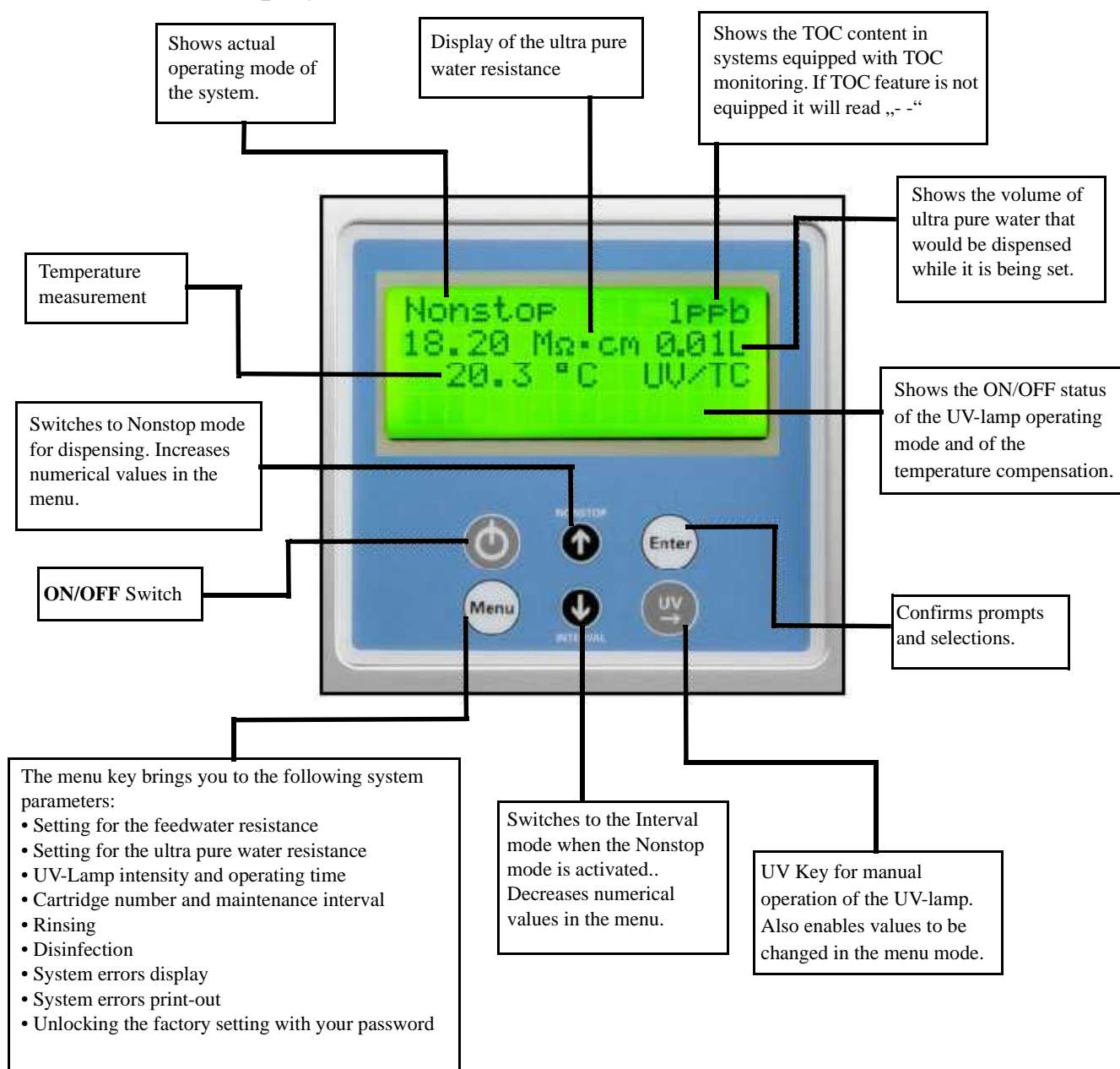
When the system successfully produces the ultrapure water quality that you require in “Nonstop” mode, press this button to return the system to the “Interval” mode.

Operating elements

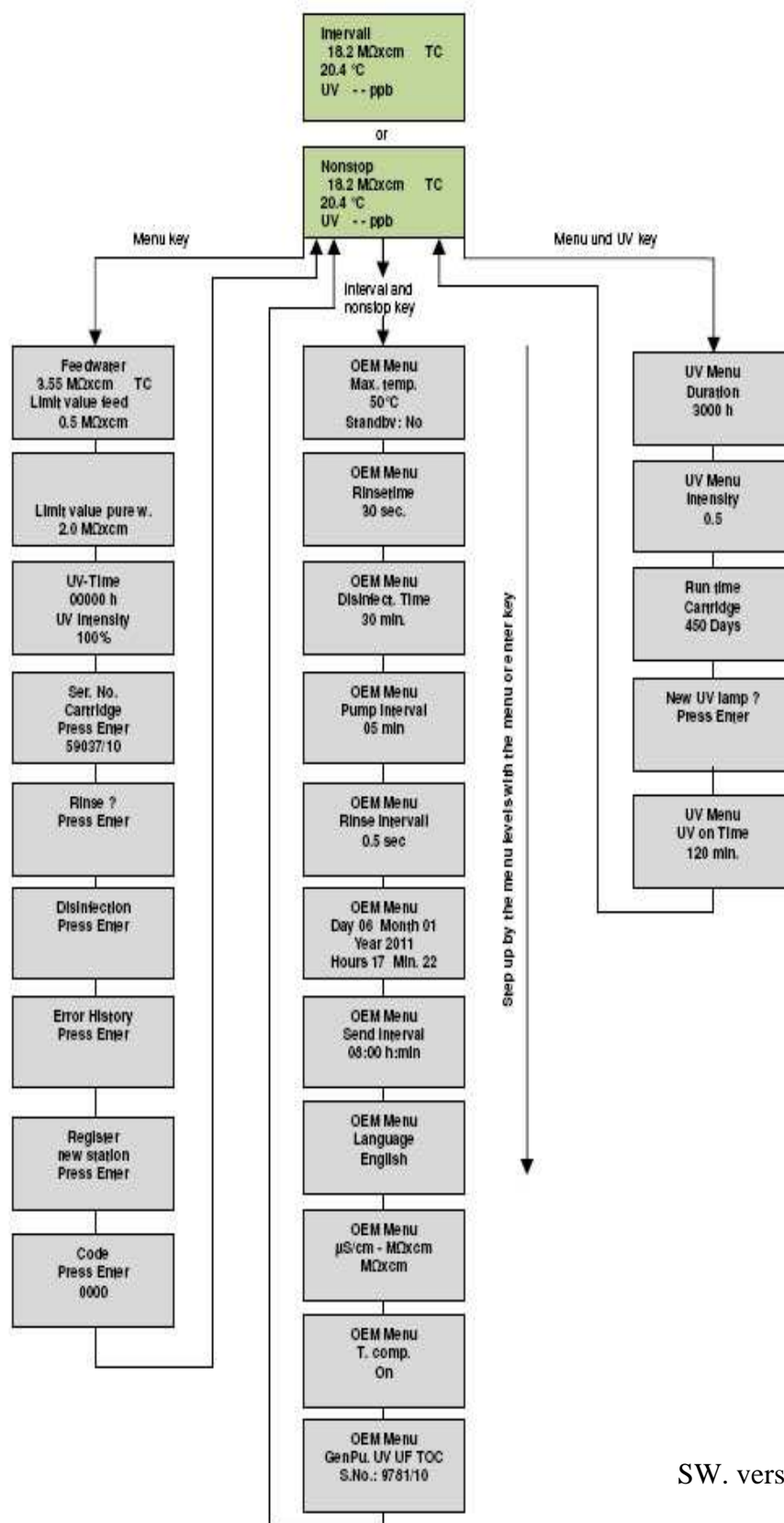
Contents

- “Description of Display” on page 55
- “Flow chart of menu system control” on page 56

Description of Display



Flow chart of menu system control



SW. version 0.0.0

11 Operating elements

Flow chart of menu system control

The system control

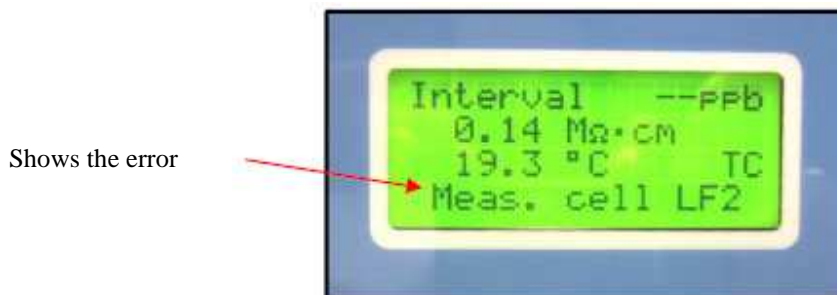
Contents

- “General information” on page 59
- “Operating modes” on page 59
- “User menu” on page 62
- “The OEM Menu” on page 68
- “Using volume control for water dispensing” on page 72
- “Printer output” on page 72

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 error 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 60](#)), whereby the green background of the display is switched on and remains in that work 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 there is no demand for the non-stop mode. The interval mode is designed to 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 “Intv.rinse time”. When the interval pump time has expired, the pump is switched off until the end of the standstill time. The rest 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 an 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 time” error message is brought to display when the limiting value set for this time is exceeded. TOC measurement is carried out during the time that the UV-lamp is glowing only.

The display shows:



Water dispensing via volumetric dispense

Ultrapure water systems which are equipped with the volumetric dispense option can dispense a preset volume of water.

As soon as the Nonstop-mode is selected, a litre volume is shown in line 2 of the display. This is the volume of ultrapure water which was last dispensed.

A single press on the Enter-button enables this volume value to be changed within the range from 0.01 to 65.5 litres by means of the arrow-buttons. The UV-button can be used to position the cursor at the particular number that you wish to change.

A second press on the Enter-button causes the volume of water that has been set to be dispensed. The liter volume shown in the display is the actual volume dispensed. Dispensing stops as soon as the set volume is reached.

Dispensing can be stopped at any time by a further press on the Enter-button. This enables small volumes to be dispensed by two successive presses on the Enter-button. One press starts dispensing and, when the wanted amount has been dispensed, a second press stops dispensing.

Volumetric dispense is supported in all program versions.

The display shows:



OFF mode

A second press on the ON/Off-button causes the display to switch off and all text messages on the display to be extinguished. button

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 guided to the next menu prompt. Settings are only possible when system control has been previously unlocked by entering a valid code number.

To simplify the change of settings, a press on the UV-button allows you to select a certain individual numbers in the numerical value. 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 error 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 M Ω xcm
Limiting value setting range:	0.1- 50.0 μ S/cm
Basic setting:	0.5 M Ω xcm

When a setting above 50 μ S/cm/ 0.02 Mxcm is entered for the limiting value, the limiting value is switched off and the word „Off“ appears in the display.

Press the Menu-button once then the display shows:



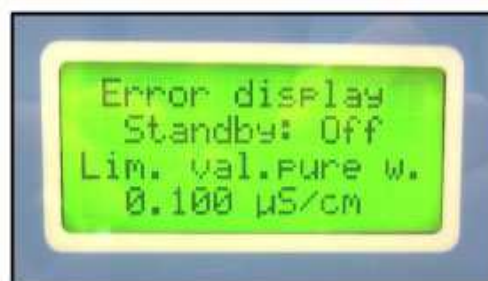
Ultrapure water limiting value

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 M Ω xcn
Limiting value setting range:	0.055- 5.000 μ S/cm
Basic setting:	10 M Ω xcn
Basic setting, fault suppression:	On

When a setting above 5.0 μ S/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:button



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:



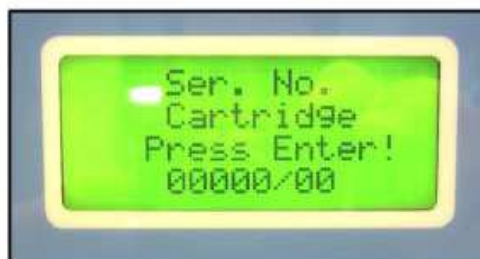
NOTE

For more details see under chapter “Change the UV-lamp” on page 80.

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:



NOTE

For more details see under chapter “Change the ultrapure cartridge” on page 71.

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:	A rectangular LCD screen with a green background and black text. The text is arranged in two lines: "Rinse?" and "Press enter". The screen is framed by a dark border.
2	Confirm rinse by putting the enter button. The rinsing is started for 30 sec.	A rectangular LCD screen with a green background and black text. The text is arranged in two lines: "Rinse" and "25 sec". The screen is framed by a dark border.

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.

During disinfection the remaining disinfection time is shown in the display.

Step	Action	Figure
1	Press the menu-button 6 times then the display shows:	
2	Confirm disinfection by putting the enter button. Change the filter cartridge with the disinfection cartridge (see under chapter “Disinfection” on page 74)	
3	Confirm with enter. The Disinfection is started for 30 min, indicating the remaining time.	

NOTE



The completely process is describe under chapter “Disinfection” on page 74” .

Error history

Confirmation of this prompt with Enter allows the fault 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 the two last recorded 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 Factory 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 13.8.

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



NOTE

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.

NOTE

Press the menu-button 10 times and you 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 programmes, 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 13.3.10).

Accessing the OEM menu.

Simultaneous presses on the INTERVAL-button and the NONSTOP-button call the OEM-menu. The display shows "OEM-Menu Press enter!". On confirming this by pressing the enter-button, the first menu point is called to be worked on. To simplify making changes, a press on the UV-button allows the position that is to be changed in a number to be selected, so that the arrow buttons can be used to replace it with any digit from 0 – 9.

A press on the menu-button takes you to the next menu point.
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 3th 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:

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 pump interval time:

In this point you can setting the pump interval time. When the system operates in the Interval mode the system is going to be recalculating every 30 min. for 5 min.

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 rinse interval 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:



Adjust the real time clock:

Basic setting: The actual date

Setting range: Month 1 - 12, Day 1 - 31, Hour 0 - 24, Minutes 0 - 60.

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



Set the sending interval:

The sending interval at which measured values and fault messages are transmitted via the RS 232 interface can be set here.

Basic setting: 1 hour

Setting range: 0.5 - 12 hours

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



Select the language:

Basic setting: German

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 $M\Omega \times cm$

Setting range: Resistance $M\Omega \times cm$, specific electrical resistance $M\Omega \text{ cm}$

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:



Using volume control for water dispensing

GenPure systems that are equipped with the option of volume control allow volume-controlled dispensing to be carried out.

As soon as the Nonstop operating mode is selected, the number of litres that were last required appears as set value in line 2 of the display.

Press once on the Enter-button if you wish to use the arrow buttons to change this set value within the permissible range of 0.01 to 65.5 litres. You can use the UV-button to position the cursor at the position where you want to change the number.

Press twice on the Enter-button if you wish to have the displayed water volume dispensed. During dispensing, the number shows the actual volume dispensed. Dispensing is stopped as soon as the set value has been reached.

Dispensing can be stopped at any time by a further press on the Enter-button.

To carry out manual dispensing of volumes smaller than the set value, first press the Enter-button twice, then press it once again when the required volume has been dispensed.

The display shows:



Printer output

By means of the printer different parameters are documented. It is differentiated between three messages.

- 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 the “Assignment Table for code numbers which unlock the system”).

Print-out:

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

Error message:

When a fault message is shown in the display, e.g. for the ultrapure water limiting value, then the fault 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 75
- “Change the ultrapure cartridge” on page 76
- “Disinfection” on page 78
- “Change the ultrafilter” on page 81
- “Structure of the UV-lamp” on page 83
- “Change the UV-lamp” on page 85
- “Change and autoclave the Final filter” on page 89

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.

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	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	12 Months	Or flow rate is noticeably slower.
Ultrafiltration membrane (only applicable for systems with a UF filter)	F3	50133980	24 Months	Or when the ultrapure water limiting value is exceeded, whichever is shorter. Longer usage can result in bacterial growth on the resin.
UV-lamp (only applicable for systems with a UV lamp)	UV1	09.2002	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

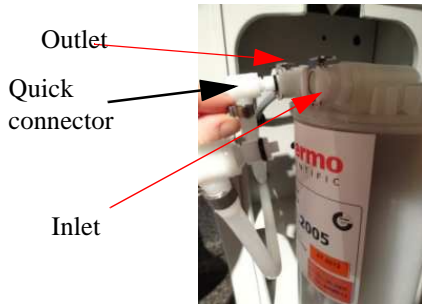


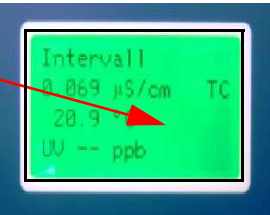





Replace the ultrapure cartridge when the maximum limiting value that you have set for the ultrapure water is exceeded or when the „New filter set“ message is shown in the display.

Step	Action	Figure
1	Switch the system off and shut off the supply of feedwater.	<p>Feedwater supply</p>
2	Remove the cartridge cover by pressing the push button.	<p>Push button</p> <p>Cartridge</p>
3	Disconnect the Quick connectors on the feedwater inlet and purified water outlet of the cartridge, close the inlet and outlet with the stoppers you have kept for later use.	<p>Outlet</p> <p>Quick connectors</p> <p>Filter cartridge</p> <p>Inlet</p>
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;">NOTE</div> <p>For sanitize your system see under chapter “Disinfection” on page 78.</p>
5	Remove the yellow stoppers from the new ultrapure cartridge and insert it into the system. Keep the yellow stoppers for the next time you have to change the cartridge.	<p>ultrapure cartridge</p>

13 Maintenance

Change the ultrapure cartridge

Step	Action	Figure
6	Plug the quick-connects correctly onto the new 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.	  
8	<div data-bbox="231 974 502 1041" style="background-color: #0056b3; color: white; padding: 5px; text-align: center; font-weight: bold;">NOTE</div> <p>For the code to perform this transaction please refer to the Code table „code lock” found in chapter „Code lock on page 67“. 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. 	<div style="display: flex; justify-content: space-around;"> <div style="text-align: center;"> <p>a.</p>  </div> <div style="text-align: center;"> <p>b.</p>  </div> </div> <div style="text-align: center; margin-top: 20px;"> <p>c.</p>  </div>
9	<div data-bbox="231 1680 502 1747" style="background-color: #0056b3; color: white; padding: 5px; text-align: center; font-weight: bold;">NOTE</div> <p>Discard at least 5 liters of water.</p>	

Disinfection



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

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

Use cleaning solutions as follows:

MICRO-Chlorine Granulate, 1 box, Catalog no. 09.2202 (Europe Emerging markets, and APAC markets)

Cleaning Solution, 1 syringe, Catalog no. CMX 25 (US and LATAM markets).



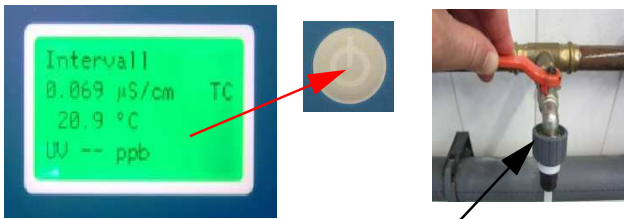
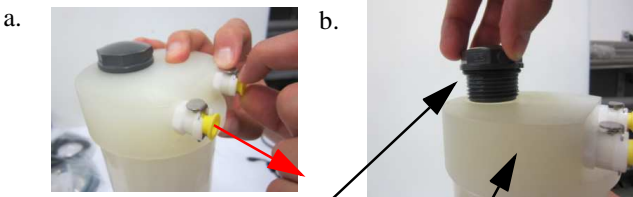
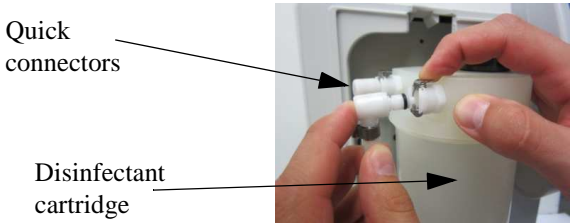
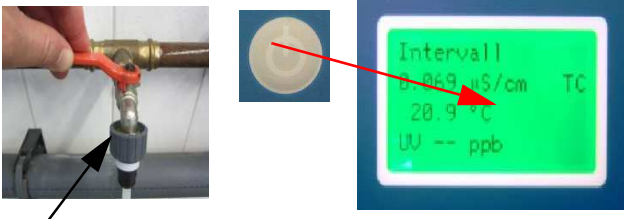
For effective disinfection the cartridge must be completely filled with distilled water.







Wear protective gloves for handling chlorine tabs or a syringe of Cleaning Solution.




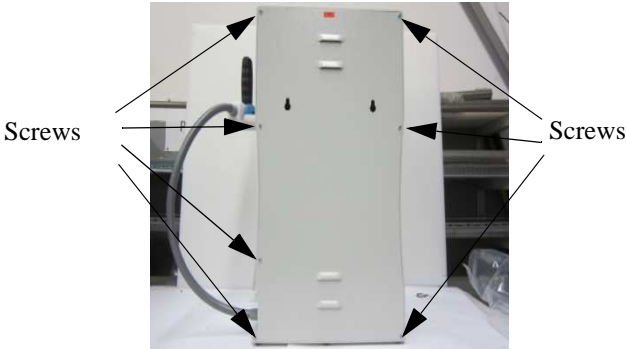
Please observe the information given in the safety data sheet supplied with Micro-Chlor disinfectant to avoid possible health hazards!


Step	Action	Figure
1	Switch the GenPure with xCAD system off and shut off the supply of feedwater. After this remove the ultrapure cartridge. NOTE See under chapter “Change the ultrapure cartridge” on page 76.	 <p>Feedwater supply</p>
2	a. Remove the yellow stoppers. b. Unscrew the stopper from the disinfectant cartridge. c. Fill the cartridge with distilled water, then empty the contents of a syringe of Cleaning solution or a can of MICRO CHLOR into the water.	 <p>a. Stopper</p> <p>b. Disinfectant cartridge</p>
3	Screw the stopper back on the disinfectant cartridge and connect the cartridge into the system. NOTE See under chapter “Change the ultrapure cartridge” on page 76 to put in the ultrapure cartridge in to the system.	 <p>Quick connectors</p> <p>Disinfectant cartridge</p>
4	Re-open the feedwater supply, switch the system on again.	 <p>Feedwater supply</p>

Step	Action	Figure
5	<p>Push the menu button until „Enter code” is displayed</p> <p>NOTE</p> <p>The Code to do this transaction please refer from the Code table under chapter „Code lock on page 67”. 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>NOTE</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>NOTE</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>NOTE</p> <p>See step 5 under chapter “Change the ultrapure cartridge” on page 76.</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>	

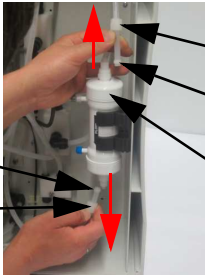
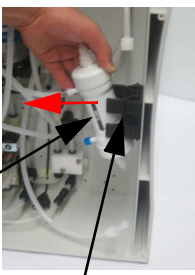



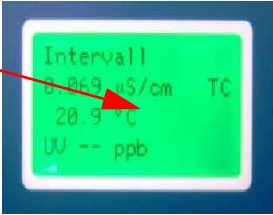
Change the ultrafilter

(applicable only for systems with UF)

Step	Action	Figure
1	Switch the GenPure with xCAD system off and shut off the supply of feedwater.	 <p>Feedwater supply</p>
2	Remove the four screws of the back panel.	

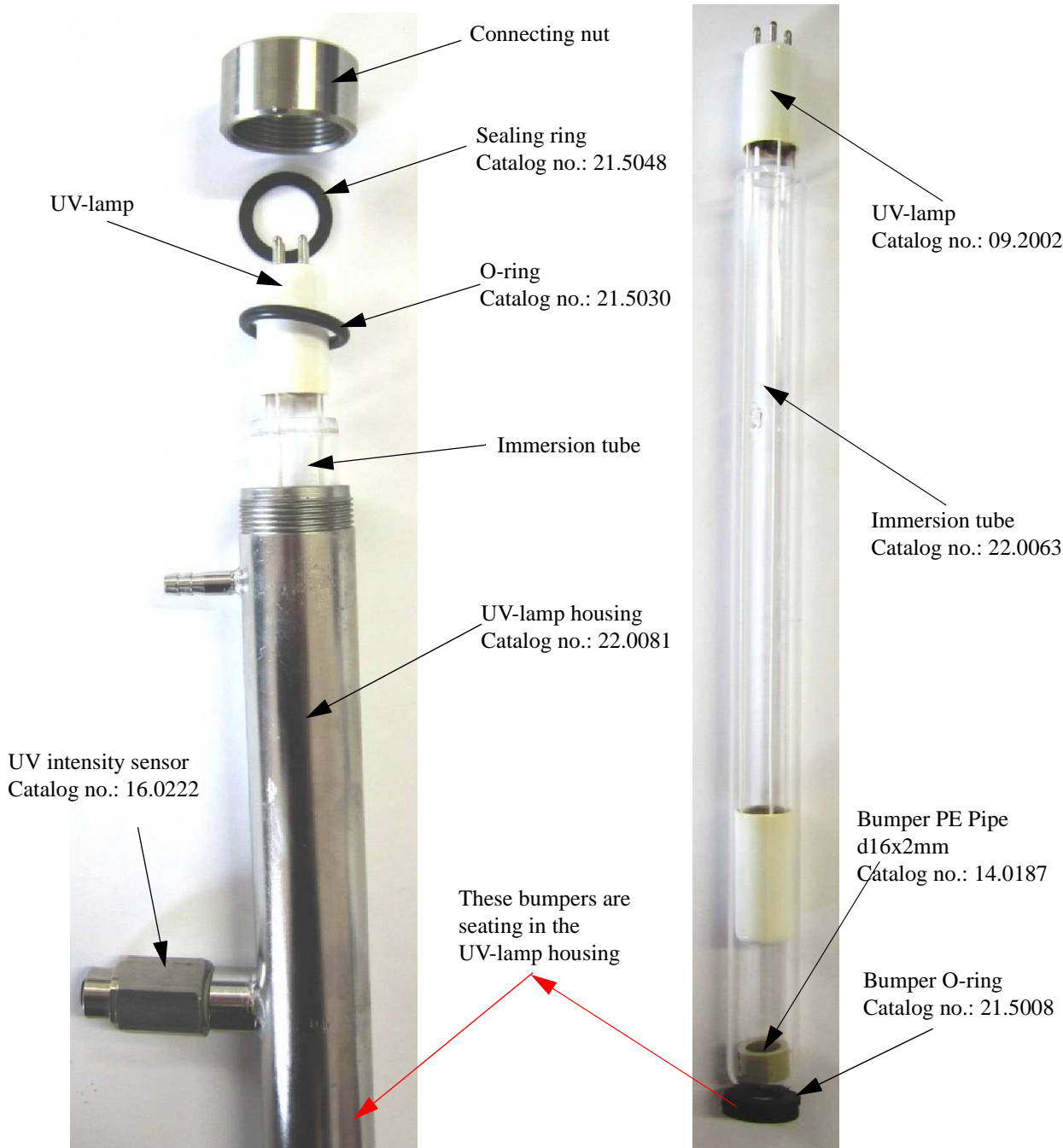


Remove carefully the back panel from the system and unscrew the yellow ground wire from the back panel.

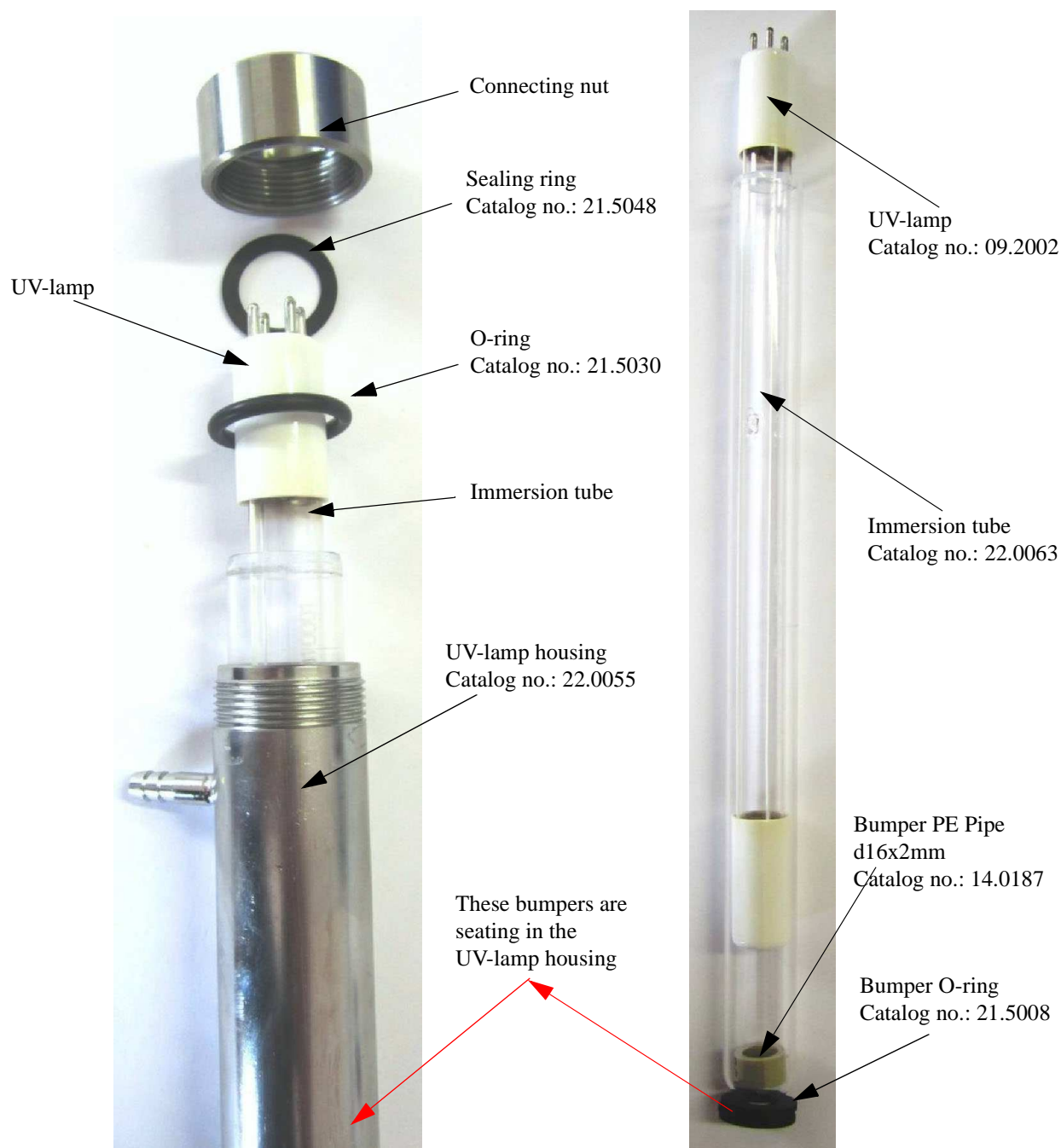
Step	Action	Figure
3	<p>a. Pull out the hoses 8 mm by unscrew the two fittings (see red arrows).</p> <p>b. After this procedure draw out the ultrafilter from the mounting clamp (see red arrow).</p> <p>c. Hold with one hand the hose and with the other hand turn in clockwise direction the ultrafilter to 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>a. </p> <p>b. </p> <p>c. </p>
	<p>NOTE</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>	
4	Reinstall the back panel, reopen the feedwater supply and switch on the system again.	<p></p> <p></p> <p></p>

Structure of the UV-lamp

UV unit with UV intensity sensor



UV unit without UV intensity sensor

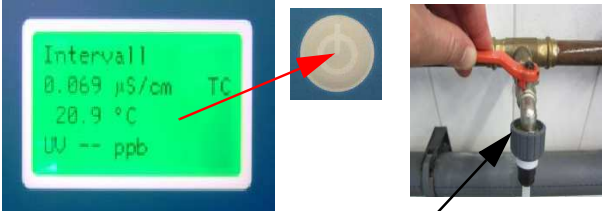
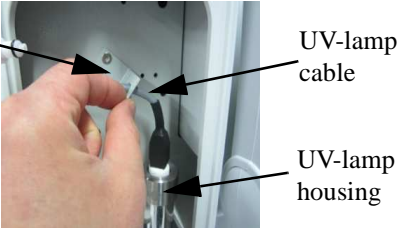
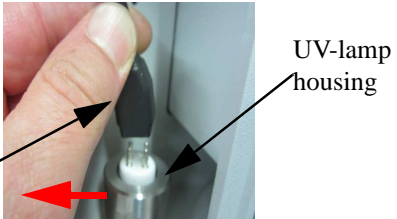


Change the UV-lamp



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

(applicable only for systems with UV lamp)

Step	Action	Figure
1	Switch the GenPure with xCAD system off and shut off the supply of feedwater.	
2	Remove the cartridge cover and take off the ultrapure cartridge.	<div><div>NOTE</div><p>See under chapter “Change the ultrapure cartridge” on page 76.</p></div>
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) and take the plug off of the UV-lamp.	

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

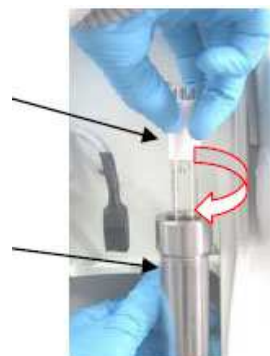
- | | |
|---|---|
| 5 | Now carefully draw the UV-lamp upwards while lightly turning it clockwise. During the replacement of a UV-lamp, great care must be taken to avoid touching the glass of the UV-lamp with fingers, to avoid dirtying of the lamp which would impair the functioning of it. |
|---|---|

NOTE

We therefore recommend that clean gloves be worn.

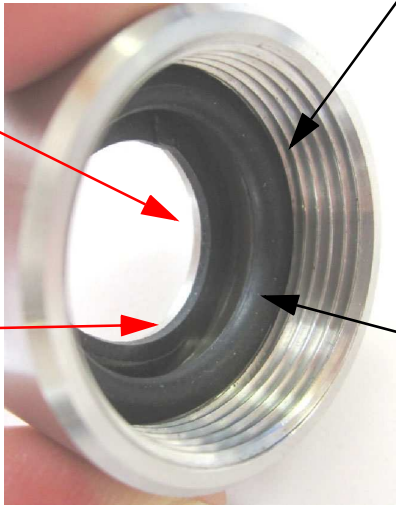
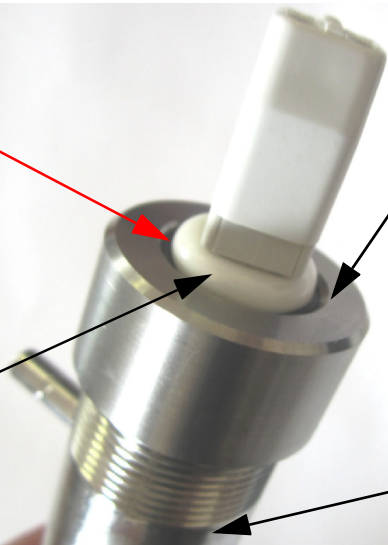
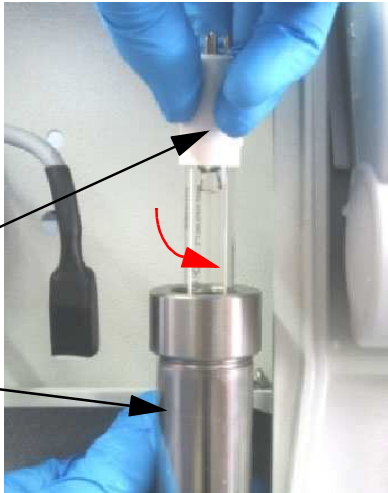
UV-lamp

UV-lamp housing



NOTE

See chapter “[Structure of the UV-lamp](#)” on [page 83](#) where is seating the sealing ring to not damage it.

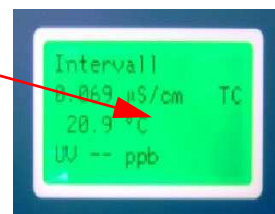
Step	Action	Figure
6	<div data-bbox="225 304 517 383" data-label="Image"> </div> <p>Ensure that the position of the sealing ring (flat o-ring at the top of the connection nut) is correct as you put in the new UV-lamp, otherwise you will have a leak. The sealing ring must be seat in the rabbet of the connecting nut (see picture a and b).</p> <p>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 remove the mounting plate.</p>	<p>a.</p>  <p>Rabbet</p> <p>Sealing ring</p> <p>Connection nut</p> <p>O-Ring</p> <p>b.</p>  <p>Sealing ring</p> <p>Connection nut</p> <p>UV lamp</p> <p>UV-lamp housing</p> <p>c.</p>  <p>UV lamp</p> <p>UV-lamp housing</p>

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

- 7 Put the cartridge cover back on (see under chapter “[Change the ultrapure cartridge](#)” on [page 76](#)), re-open the feed water supply and switch the system on again.



Feedwater supply



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

NOTE

The Code to do this transaction please refer to the Code table under chapter „[Code lock](#) on [page 67](#)”. You need a level 3 code.

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

a.



b.



c.

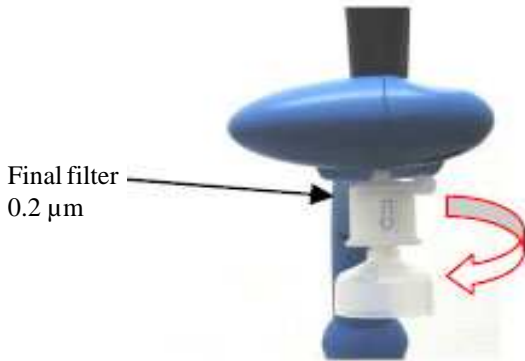
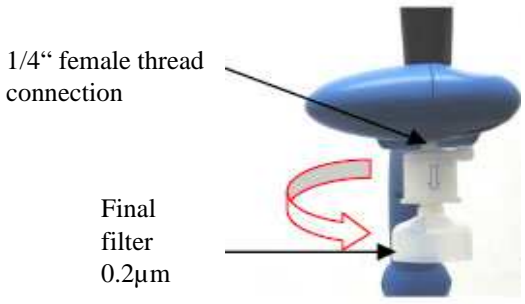


NOTE

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.

Change and autoclave the Final filter

Step	Action	Figure
1	Unscrew the blocked or used final filter by turning it in clockwise direction.	
2	Unpacking the new Final filter and screw in the it in the dispensing valve outlet (R 1/4" female thread).	

Autoclave the Final filter

NOTE

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 10 times. When the sterilization is finished screw in the final filter back in the 1/4" female thread connection (see chapter [“Change and autoclave the Final filter”](#) on [page 89](#)).

NOTE

If you 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 94](#) or change with a new one.

13 Maintenance

Change and autoclave the Final filter

Waste disposal

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

NOTE

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

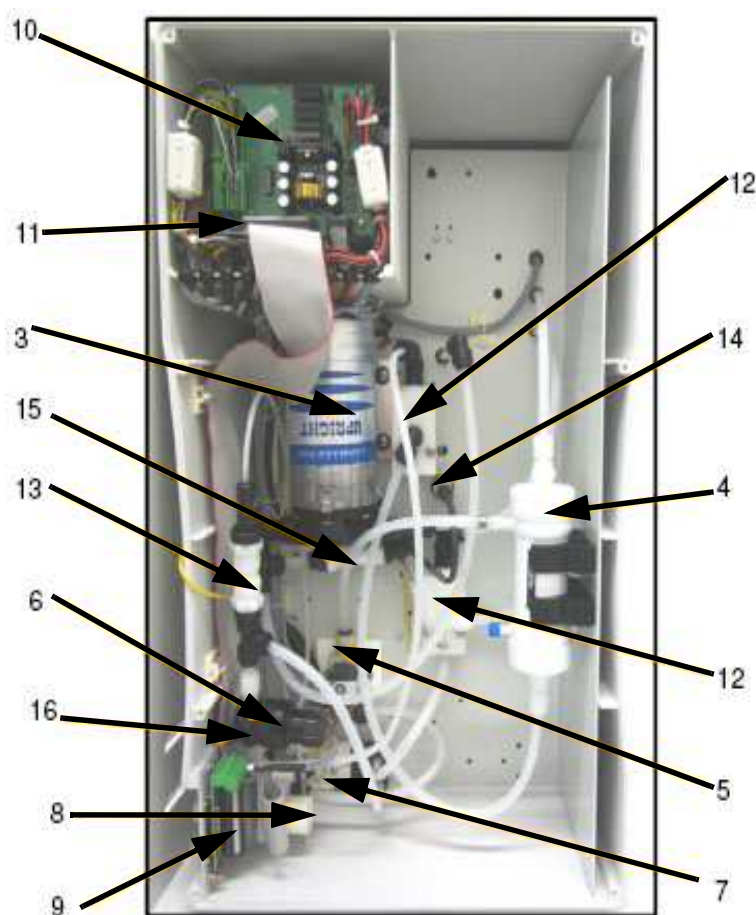
Error	Cause	Remedy
The system does not start	<ul style="list-style-type: none"> No supply of power 	<ul style="list-style-type: none"> Provide power
Dispensing not possible	<ul style="list-style-type: none"> Feedwater tap is closed Feedwater and rinse water connections are mixed up Feedwater pressure < 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 MΩxcm	<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 ultrapure 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
Dispensed amount is too small	<ul style="list-style-type: none"> UF-Module blocked Pre-pressure too low Internal pressure too low Volumetric Dispense out of Tolerance 	<ul style="list-style-type: none"> Replace UF-module Increase the pre-pressure Readjust pressure reducer Contact Thermo for volume calibration

Error	Cause	Remedy
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: „Limit value feed“	<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: „Lim. va.pure w.“	<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: „UV-time“	<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: „UV-intensity“	<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 with a new UV-lamp Clean the UV-sensor Check and set the limiting value
Error message: „max. Temperature“	<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: „Measuring cell LF1“	<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 MΩxcm” on page 94
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 95

Error	Cause	Remedy
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 and consumables

GenPure



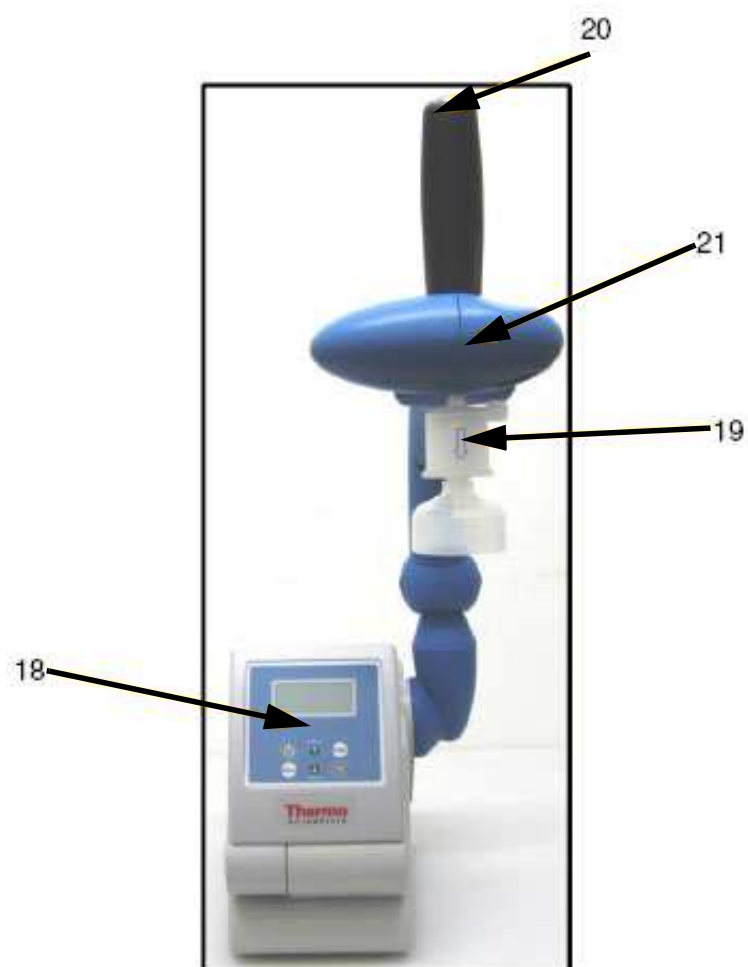
Parts marked with an „x“ are wear parts (exchange is performed by service people)

No.	Flow chart no.	Designation	Catalog no.
1	F1	Ultrapure cartridge	09.2005
2	UV1	Replacement UV lamp	09.2002
3	P1	Pressure booster pump	19.0050 x
4	F3	Ultrafiltration module (optional)	50133980
5	V4	Rinsing solenoid valve	50131190 x
6	V1	Pressure reducer	50133985
7	FIS400	Flowmeter	15.0100
8	QIA300	Conductivity measuring cell, feedwater	16.0126
9		Adapter board	16.0364
10		Microprocessor-system control, interface	50131346
11		Booster for UV lamp (optional)	22.0088
12	QIA303 TIA500	Conductivity measuring cell, ultra pure water Conductivity measuring cell, temperature sensor	50133992
13	V2	Check valve	15.0009
14	QIA301	UV-Intensity sensor (optional)	16.0222
15	V5	Check valve	15.0019
16		Fuseholder for glas tube fuse, 5x20mm Glas tube fuse, 5x20mm, 3.15 A, slow	50133979 50131758
17		Table top power pack, 24V DC (not shown)	50134196

NOTE

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.

xCAD



Parts marked with an „x“ are wear parts (exchange is performed by service people)

No.	Flow chart no.	Designation	Catalog no.
18		Microprocessor-system control	26.0025
19	F2	Final filter 0.2 µm	09.1003 x
20		Press button	16.0370
21	V3	Solenoid valve	15.0101 x
22		Extension cable SUB-D, 25 pin, GenPure/xCAD (not shown)	16.0375

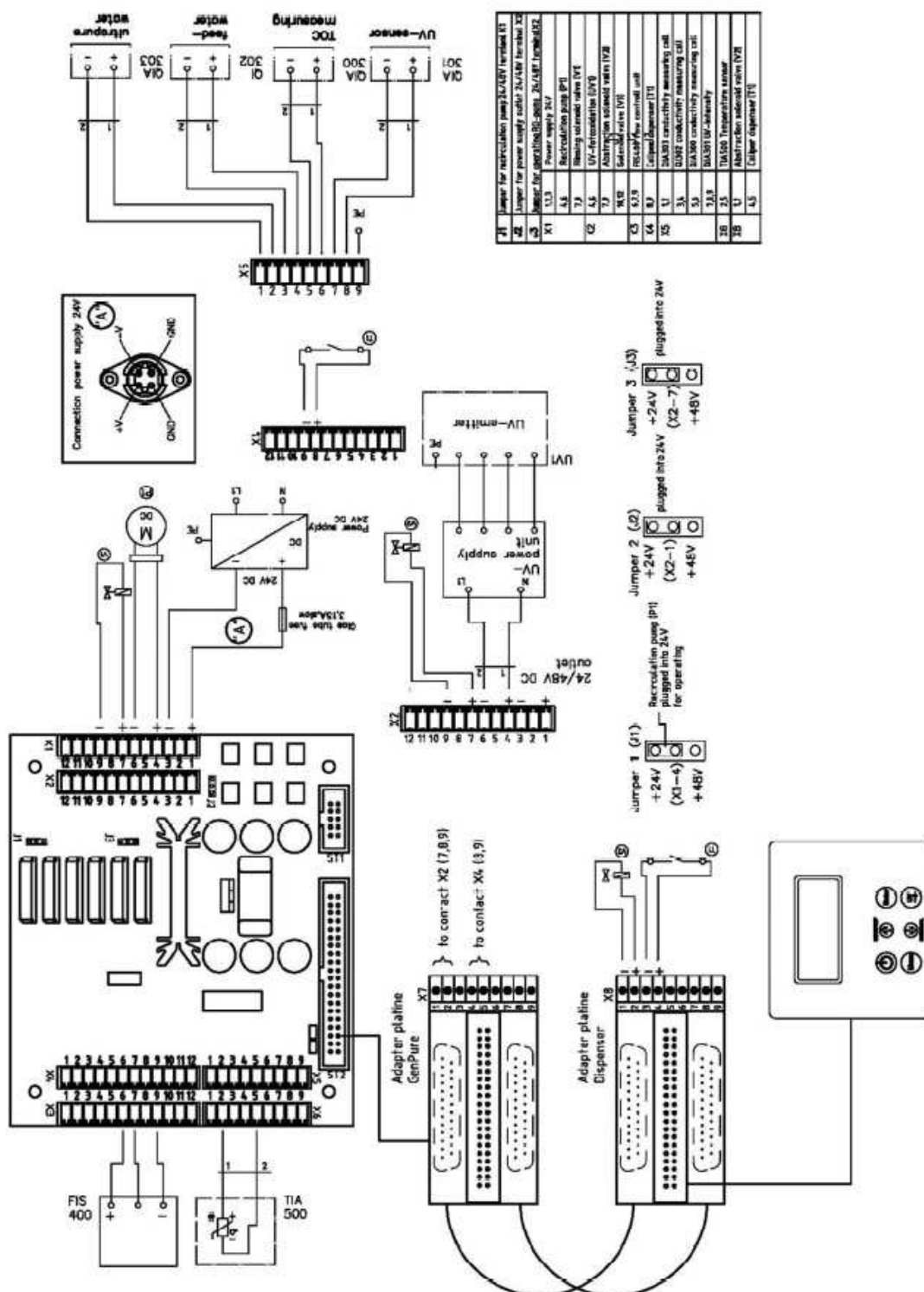
Consumable materials

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
Disinfection agent, MICRO-Chlor (pack of 12 cans, Europe only)	09.2202
Cleaning Solution, 1 syringe (US-market only)	CMX25
Printer	09.2207
Ion exchanger DI 1500	02.1500
DI 1500 hose kit for new installations	04.1690

Terminal assignment



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 ultrapure 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

© 2013 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 [102](#)
 Autoclave
 Final filter [90](#)
 Available GenPure Systems [12](#)

C

Change
 Final filter [89](#)
 Ultrafilter [81](#)
 UV-lamp [85](#)
 Code message [78](#)
 Complaints [6](#)
 Consumable materials [102](#)
 Contact information [110](#)

D

Data transmission via the RS 232 interface [77](#)
 Dimension and weight [22](#)
 Disinfection [78](#)

E

Examination on receipt [6](#)
 Explanatory notes [v](#)
 Extent of assembly kit [110](#)
 Extent of delivery [19](#)

F

Filter cartridge
 Change [76](#)
 Final
 Autoclave [90](#)
 Final filter
 Change [90](#)
 Flow chart control unit [56](#)
 Flow charts [42](#)
 GenPure Standard [43](#)
 GenPure UF [44](#)
 GenPure UV [45](#)
 GenPure UV/UF [46](#)
 GenPure UV-TOC [47](#)

GenPure UV-TOC/UF [48](#)

I

Information, legal [iv](#)
 Installation [25](#)
 Installation area [23](#)
 Intended use [115](#)

L

Legal Information [iv](#)

M

Maintenance [74](#)
 Intervals [75](#)
 Records [108](#)
 Mounting the power pack [39](#)

O

OEM menu [68](#)
 Changing the disinfection time [69](#)
 Language selection [71](#)
 Rinsing time [68](#)
 Setting the interval pump time [69](#)
 Setting the sending interval [70](#)
 Switching temperature compensation off [71](#)
 Switching units [71](#)
 Operating elements [54](#)
 Operating modes [59](#)
 Interval operating mode after switching on [59](#)
 Interval operation [60](#)
 Non-stop mode [60](#)
 OFF mode [62](#)
 UV-Lamp [60](#)
 Water dispensing via volume control [61](#)

P

Packaging [6](#)
 Packing for return shipment [6](#)
 Printer output [72](#)

Index:

Print-out [71](#)
Putting into operation [52](#)

R

Replacement parts [98](#)
Return shipment [6](#)

S

Safety precautions [7](#)
Setting
 Interval pump time [69](#)
 Sending interval [70](#)
Specifications [17](#)
Standard message [72](#)
Start your system into operation [52](#)
System
 Functions [54](#)
 Putting into operation [52](#)
System control [63](#)
 General information [59](#)

T

Technical specifications [17](#)
 Airborne sound emission [17](#)
 Ambient conditions [17](#)
 Cell constants of the measuring cells [68](#)
 Connectors for water [17](#)
 Demands the feedwater must fulfill [17](#)
 Electrical connections / external switched mode power supply [17](#)
 Materials of parts which contact water [17](#)
 Product water quality [17](#)
Terminal assignment [1106](#)
Transport [5](#)
Trouble shooting [94](#)

U

Ultrafilter
 Change [81](#)
User menu [62](#)
 Entering a code number [67](#)
 Error history [66](#)
 Feedwater measured value and limiting value [62](#)
 Ultrapure cartridge serial number [76](#)
 Print out of Data [72](#)
 Rinsing the ultrafilter [64](#)
 Ultrapure water limiting value [63](#)
 UV Intensity and operating time [63](#)
UV-lamp
 Change [85](#)
 Structure [83](#)

W

Warranty [iv](#)
Waste disposal [92](#)



Thermo Fisher Scientific,
San Jose, CA USA is ISO Certified.

thermoscientific.com

© 2013 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
S C I E N T I F I C
Part of Thermo Fisher Scientific