

Operating Instructions

Original Operating Instructions

Arium® Mini Plus - Arium® Mini Plus | Extend

Arium® Mini

Arium® Mini Essential - Arium® Mini Essential | Extend

H2O-MA-... | H2O-MAR-... | H2O-MM-... | H2O-MU-... | H2O-MUR-...

Ultrapure Water System



1000139721



SARTORIUS

Contents

1 About this Document	5	5 Installation	30
1.1 Validity	5	5.1 Equipment Supplied	30
1.2 Symbols Used	5	5.2 Prerequisites for Installation at the Installation Location	31
1.2.1 Warnings in Action Descriptions	5	5.3 Unpacking and Setting up the Device	32
1.2.2 Other Symbols	5	5.3.1 Removing the Left Side Cover	32
2 Safety Notes	6	5.3.2 Removing Parts	33
2.1 Intended Use	6	5.4 Mounting the Dispensing Unit	33
2.2 Personnel Qualification	6	5.4.1 Connecting the Tube Unit	34
2.3 Significance of These Instructions	7	6 Startup	35
2.4 Functionality of the Device	7	6.1 Connecting the AC Adapter	35
2.5 Electrical Equipment	8	6.2 Configuring Device Settings (Startup)	36
2.5.1 Damage to the Electrical Equipment of the Device	8	6.2.1 Setting the Language	36
2.5.2 Working on the Electrical Equipment of the Device	8	6.2.2 Launching Startup Mode	37
2.5.3 Power Supply and Power Cord	8	6.2.3 Setting the Date and Time	37
2.6 Accessories, Consumables and Spare Parts	8	6.2.4 Setting the Displayed Values	38
3 Device Description	9	6.2.5 Concluding Device Settings (Startup)	38
3.1 Device Overview	9	6.3 Inserting Pretreatment Cartridge (only Arium® Mini Plus)	39
3.1.1 Hand-held Part	10	6.4 Inserting Ultrapure Water Cartridge	40
3.2 Electrical Connections	11	6.5 Inserting Bag (only Arium® Mini Plus or Arium® Mini)	41
3.3 Water Connections	12	6.6 Connecting Tubing	42
3.4 Tubing	13	6.6.1 Connecting Feed Water Tubing (only Arium® Mini Plus or Arium® Mini)	42
3.4.1 Feed Water Tubing	13	6.6.2 Connecting Tank Filling Tubing (only Arium® Mini)	43
3.4.2 Tank Filling Tubing	13	6.6.3 Connecting Drain Water Tubing (only Arium® Mini Plus or Arium® Mini)	44
3.4.3 Drain Water Tubing	13	6.6.4 Connecting the Tank Outlet Tubing (only Arium® Mini Plus or Arium® Mini)	45
3.4.4 Tank Outlet Tubing with Ball Cock	13	6.7 Rinsing Pretreatment Cartridge (only Arium® Mini Plus)	45
3.4.5 Dispense Tube	13	6.8 Rinsing the Ultrapure Water Cartridge	46
3.5 Ultrapure Water Treatment	14	6.8.1 Filling of the Bag (only Arium® Mini Plus and Arium® Mini)	46
3.5.1 System Setup	14	6.8.2 Performing a Rinse Cycle	47
3.5.2 Feed Water Intake	16	6.8.3 Re-filling the Bag During the Rinse Cycle (only Arium® Mini)	48
3.5.3 Treatment Stages	16	6.8.4 Completing the Rinsing Process	49
3.5.4 Final Purification Stage	17	6.9 Connecting the Final Filter	49
3.5.5 Circulation of the Ultrapure Water	17	6.10 Rinsing the Final Filter	50
4 Operating Design	18		
4.1 Menu	18		
4.2 Dispensing Mode	19		
4.3 Message Display	20		
4.4 Messages in Dispensing Mode	21		
4.5 Message List	22		
4.6 Numeric Keypad	23		
4.7 Volume Selection (Only for Devices With a Dispensing Unit)	24		
4.8 Navigating the Menus	25		
4.9 Menu Structure	26		
4.10 Parameters of the "Settings" Menu	28		

7 Operation	51	9 Malfunctions	73
7.1 Switching the Device On and Off	51	9.1 Error Messages	73
7.2 Dispensing Ultrapure Water	51	9.2 Warning Messages	76
7.2.1 Preparing to Dispense Ultrapure Water	51	9.3 Additional Faults	78
7.2.2 Dispensing Ultrapure Water Manually	52	10 Storage and Shipping	79
7.2.3 Dispensing Ultrapure Water with Volume-control	53	10.1 Storage	79
7.3 Dispensing Ultrapure Water From the Bag (Only With Arium® Mini Plus or Arium® Mini)	55	10.2 Returning Device and Parts	79
7.4 Opening the Menu	55	11 Decommissioning	80
7.5 Manually Fill the Bag (Only for Arium® Mini)	56	12 Transport	80
7.5.1 Starting Manual Filling via the Wizard	56	12.1 Transporting the Device	80
7.5.2 Starting Manual Filling in Menu	56	13 Disposal	81
7.6 Activating or Deactivating Standby Mode	58	13.1 Information on Decontamination	81
7.6.1 Automatic ECO Mode	58	13.2 Disposing of Device and Parts	81
7.7 Changing System Settings	58	13.2.1 Information on Disposal	81
7.8 Performing a Volume Adjustment	59	13.2.2 Disposal	82
8 Cleaning and Maintenance	60	14 Technical Specifications	83
8.1 Cleaning	60	14.1 Power Supply	83
8.1.1 Cleaning the Display	60	14.2 Ambient Conditions	83
8.1.2 Cleaning the Device Housing	60	14.3 Safety of Electrical Equipment	84
8.2 Maintenance Schedule	61	14.4 Electromagnetic Compatibility	84
8.3 Opening the "Care" Menu	61	14.5 Product Water Quality	85
8.4 Displaying Reminders	62	14.5.1 Arium® Mini Plus and Arium® Mini Plus Extend	85
8.5 Replacing Consumables	62	14.5.2 Arium® Mini, Arium® Mini Essential and Arium® Mini Essential Extend	86
8.5.1 Selecting Consumables for Replacement	62	14.6 Feed Water Quality	86
8.5.2 Replacing Bag (only Arium® Mini Plus or Arium® Mini)	63	14.6.1 Arium® Mini Plus and Arium® Mini Plus Extend	86
8.5.3 Replacing Pretreatment Cartridge (only Arium® Mini Plus)	64	14.6.2 Arium® Mini	87
8.5.4 Replacing Ultrapure Water Cartridge	65	14.6.3 Arium® Mini Essential and Arium® Mini Essential Extend	87
8.5.5 Replacing the UV Lamp	66	14.7 Device Properties	88
8.5.6 Changing Final Filter	69	15 Consumables	89
8.6 Enabling, Disabling or Configuring Reminders for Replacing Final Filters	70	16 Sartorius Service	89
8.7 Carrying Out Depressurization	71	17 Conformity	89
8.8 Carrying Out Venting	71	17.1 EC/EU Declaration of Conformity	89
8.9 Performing the Final Rinsing Procedure	72	17.2 UK Declaration of Conformity	89

1 About this Document

1.1 Validity

These instructions are part of the device. These instructions apply to the device in the following versions:

Device	Model name
Devices without a dispensing unit	
Arium® Mini	H2O-MM-UV-T H2O-MM-T H2O-MM-UV-T-US H2O-MM-T-US
Arium® Mini Plus	H2O-MA-UV-T H2O-MA-UV-T-US H2O-MA-T H2O-MA-T-US
Arium® Mini Essential	H2O-MU-UV-T H2O-MU-T H2O-MU-UV-T-US H2O-MU-T-US
Devices with a dispensing unit	
Arium® Mini Plus Extend	H2O-MAR-UV-T H2O-MAR-T
Arium® Mini Essential Extend	H2O-MUR-UV-T H2O-MUR-T

1.2 Symbols Used

1.2.1 Warnings in Action Descriptions

WARNING

Denotes a hazard that may result in death or serious injury if it is **not** avoided.



CAUTION

Denotes a hazard that may result in moderate or minor injury if not avoided.

NOTICE

Denotes a hazard that may result in property damage if it is not avoided.

1.2.2 Other Symbols

-  Required action: Describes activities that must be performed.
-  Result: Describes the result of the activities carried out.
- [] Text inside brackets refers to control and display elements.

Figures on the Operating Display

Depending on the device configuration, the figures depicting the device and operating display may differ slightly from the supplied device. The variants shown in these instructions are examples.

2 Safety Notes

2.1 Intended Use

The device is a water treatment system, which produces ultrapure water of stage "ASTM type 1" for a daily requirement of up to 10 liters. In the case of devices with a dispensing unit, the dispensing of the ultrapure water is also controlled via the hand-held part of the dispensing unit.

The device is intended exclusively for use in accordance with these instructions. Any further use beyond this is considered **improper**.

If the device is **not** used properly: The protective systems of the device may be impaired. This can lead to unforeseeable personal injury or property damage.

Operating Conditions for the Device

Do **not** use this device in potentially explosive environments. The device may only be used indoors.

The device may only be used with the equipment and under the operating conditions described in the Technical Data of these instructions.

Modifications of the Equipment

You may **not** modify the device or make any technical changes on your own. Any retrofitting or technical changes to the device are only permitted with prior written permission by Sartorius.

Foreseeable Misuse

Using the device is only safe when it is operated in accordance with its intended use. The following applications, for example, are **not** permitted:

- Operation outside the permissible ambient conditions (see Chapter "14.2 Ambient Conditions," page 83) e.g. extreme temperatures, chemical vapors, moisture, shock, vibration or strong electromagnetic fields
- Carrying out unauthorized modifications or other technical changes on the device
- Connecting unsuitable devices
- Installation of unauthorized items on or in the device

2.2 Personnel Qualification

These instructions are addressed to the target groups mentioned below. All persons working on the device must possess the stated knowledge and authorizations.

If **no** qualifications are indicated for the actions described in these instructions: The actions described are addressed to the "user" target group.

If individual actions must be carried out by other target groups or by Sartorius Service personnel: The qualification required will be indicated in the description of the action.

Target group	Knowledge / responsibilities
User	<p>The user is familiar with the operation of the device and the associated work processes. He knows the dangers that can occur when working with the device and can avoid these dangers.</p> <p>The user has been trained in the operation of the device.</p> <p>The training is carried out by the operating engineer/laboratory manager or the operator of the device.</p>
Operating engineer/ laboratory manager	<p>The operating engineer/laboratory manager makes decisions about the use and parameterization of the device.</p> <p>The operating engineer/laboratory manager is trained in the operation of the device.</p> <p>The training is performed by Sartorius Service or the operator.</p>
Operator	<p>The operator of the device is responsible for compliance with safety requirements and workplace safety regulations.</p> <p>The operator must ensure that all persons who work with the device have access to the relevant information and have been instructed in work with the device.</p>

2.3 Significance of These Instructions

Failure to follow these instructions can have serious consequences, e.g. exposure of individuals to electrical, mechanical or chemical hazards.

- ▶ Before working with the device: Read the instructions carefully and completely.
- ▶ If the instructions are lost: Request a replacement or download the latest instructions from the Sartorius website (www.sartorius.com).
- ▶ The information contained in these instructions must be made available to all individuals working with the device.

2.4 Functionality of the Device

A damaged device or worn parts may lead to malfunctions or cause hazards which are difficult to recognize.

- ▶ Only operate the device when it is in proper working order.
- ▶ Comply with the maintenance intervals (for intervals and maintenance tasks, see Chapter “8.2 Maintenance Schedule,” page 61).
- ▶ Have any damage repaired immediately by Sartorius Service.

2.5 Electrical Equipment

2.5.1 Damage to the Electrical Equipment of the Device

Damage to the electrical equipment of the device, e.g. damage to the insulation, can be life-threatening. Contact with parts under voltage represents a direct danger to life.

- ▶ If the electrical equipment of the device is defective, immediately switch off the power supply and contact Sartorius Service.
- ▶ Keep live parts away from moisture. Moisture can cause short circuits.
- ▶ Make sure that the power connection is equipped with a ground lead.

2.5.2 Working on the Electrical Equipment of the Device

Any work on or modifications to the electrical equipment of the device may only be carried out by Sartorius Service personnel. The device may only be opened by Sartorius Service personnel.

2.5.3 Power Supply and Power Cord

Serious injury can result, e.g. from electric shocks, if an unsuitable power supply or an unsuitable / inadequately dimensioned power cords is used.

- ▶ Use only original power supplies and original power cords.
- ▶ **Do not** use inadequately dimensioned power cords.
- ▶ If the power supply or the power cord must be replaced:
Contact Sartorius Service. **Do not** repair or modify the power supply or power cord.

2.6 Accessories, Consumables and Spare Parts

The use of unsuitable accessories, consumables and spare parts can affect the functionality of the device, be hazardous and have the following consequences:

- Risk of injury to persons
 - Damage to the device
 - Malfunctions of the device
 - Total failure of the device
- ▶ Only use accessories, consumables and spare parts from Sartorius. Sartorius can provide information on operational quality upon request.
 - ▶ Only use accessories, consumables and spare parts that are in technically perfect condition.

3 Device Description

3.1 Device Overview

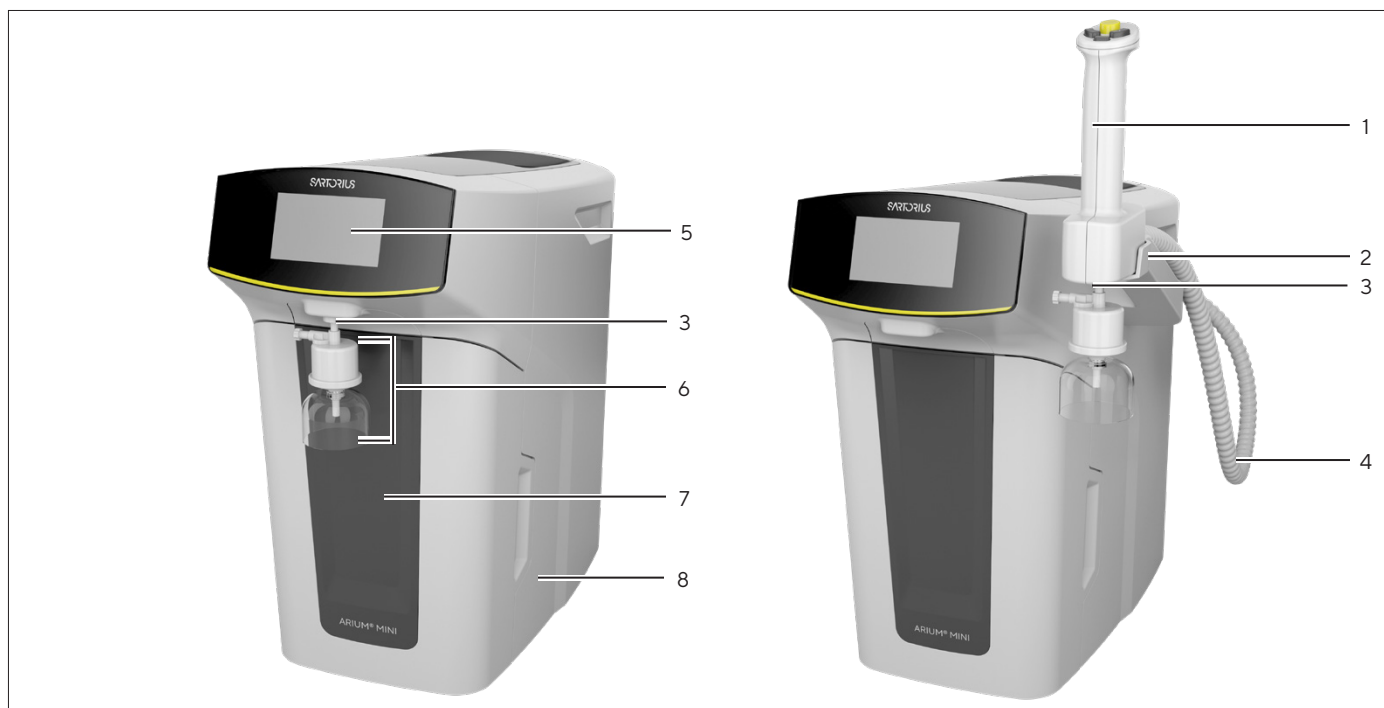


Fig.1: Arium® Mini Plus and Arium® Mini Plus | Extend (example)

No.	Description
1	Hand-held Part
2	Bracket
3	Water outlet
4	Connection tube
5	Display with touch function
6	Final filter
7	Front cover
8	Side cover

3.1.1 Hand-held Part

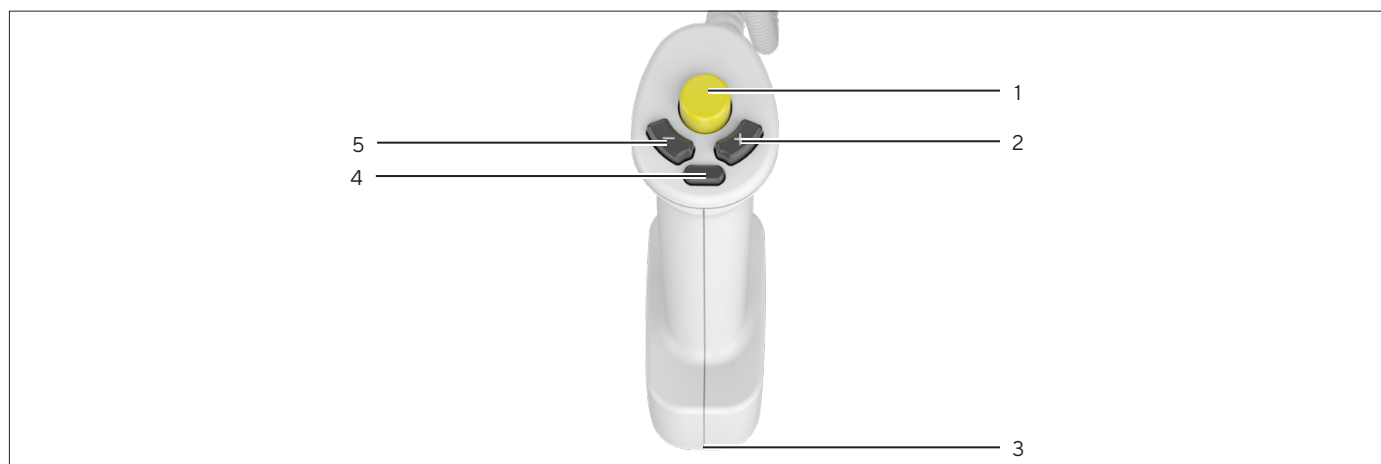


Fig.2: Handteil der Dosiereinheit (Beispiel)

Pos.	Designation	Explanation
1	[Start Stop] button	Starts or stops the dispensing.
2	[Plus] button	Increases the amount of water to be dispensed by a previously specified value.
3	Water outlet	Used for the connection of the final filter.
4	[Volume] button	Starts the selection of the water amount for volume-controlled dispensing.
5	[Minus] button	Decreases the amount of water to be dispensed by a previously specified value.

3.2 Electrical Connections

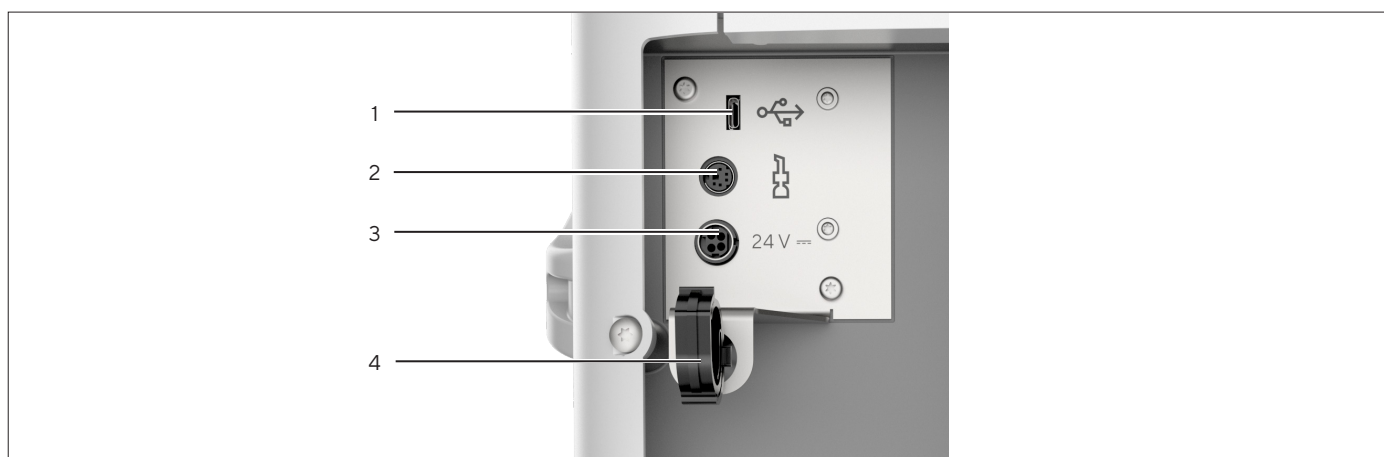


Fig.3: Electrical connections on the Arium® Mini Plus (example)

No.	Description	Explanation
1	“USB” connection	For Sartorius Service
2	Dispensing unit connection	For the electrical connection of the dispensing unit.
3	“Power” connection	For connection of the AC adapter (power supply)
4	Clip	Only for devices with a dispensing unit: Secures the tube connection to the device.

3.3 Water Connections

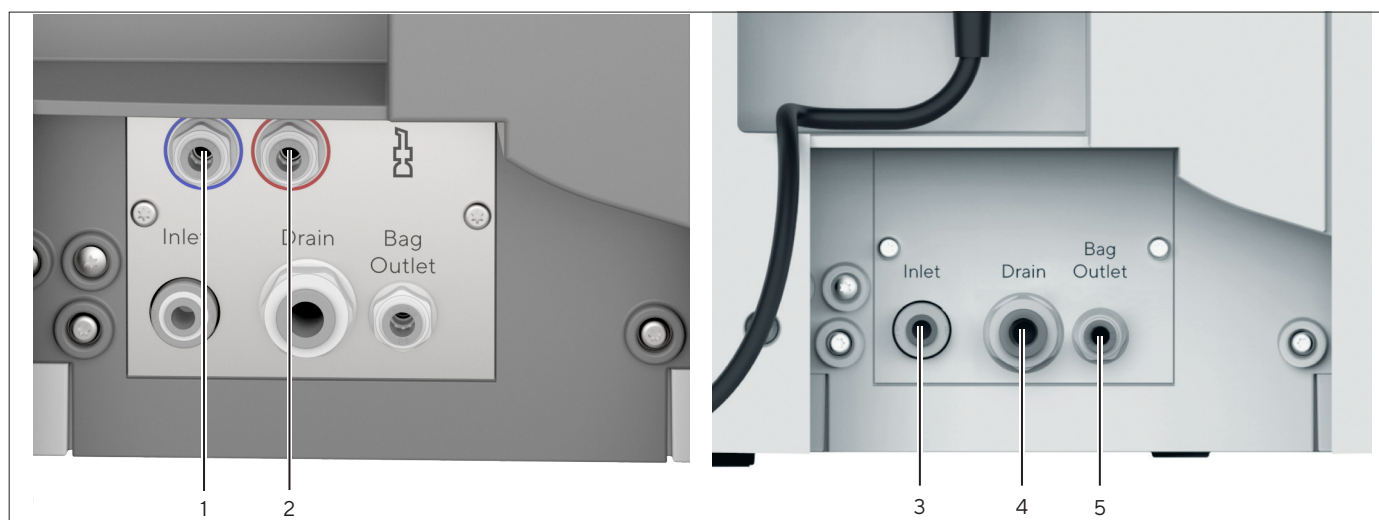


Fig.4: Water connections on the Arium® Mini Plus | Extend and Arium® Mini Plus (example)

Pos.	Designation	Explanation
1	“Blue” dispensing unit connection	For devices with a dispensing unit: For connecting the feed water tube to the hand-held part.
2	“Red” dispensing unit connection	For devices with a dispensing unit: For connecting the return water tube from the hand-held part.
3	“Inlet” connection	For the connection of the tank filling tube or inlet water tube.
4	“Drain” connection	Only Arium® Mini Plus with or without dispensing unit and Arium® Mini: For the connection of the waste water tube.
5	“Bag outlet” connection	Only Arium® Mini Plus with or without dispensing unit and Arium® Mini: For the connection of the tank outlet tube.

3.4 Tubing

3.4.1 Feed Water Tubing



The feed water tubing is used to supply feed water to the device:

- Arium® Mini Plus with or without dispensing unit: Supply of tap water
- Arium® Mini Essential with or without dispensing unit: Supply of pretreated water

The feed water tubing is marked “Inlet”.

3.4.2 Tank Filling Tubing



The tank filling tubing serves to feed pretreated water from an external tank to the device (only Arium® Mini).

The tank filling tubing is marked “Inlet”.

3.4.3 Drain Water Tubing



The drain water tubing is used to drain unpurified water and rinse water out of the device.

The drain water tubing is marked “Drain”.

3.4.4 Tank Outlet Tubing with Ball Cock



Pure water can be dispensed without pressure directly from the bag via the tank outlet tubing with ball cock. This option can be used to empty the bag when performing maintenance or to dispense pretreated water directly from the bag for further use.

The tank outlet tubing is marked “Bag Outlet”.

3.4.5 Dispense Tube



The dispense tube is used for rinsing functions during startup and maintenance as well as when dispensing larger volumes of water.

3.5 Ultrapure Water Treatment

3.5.1 System Setup

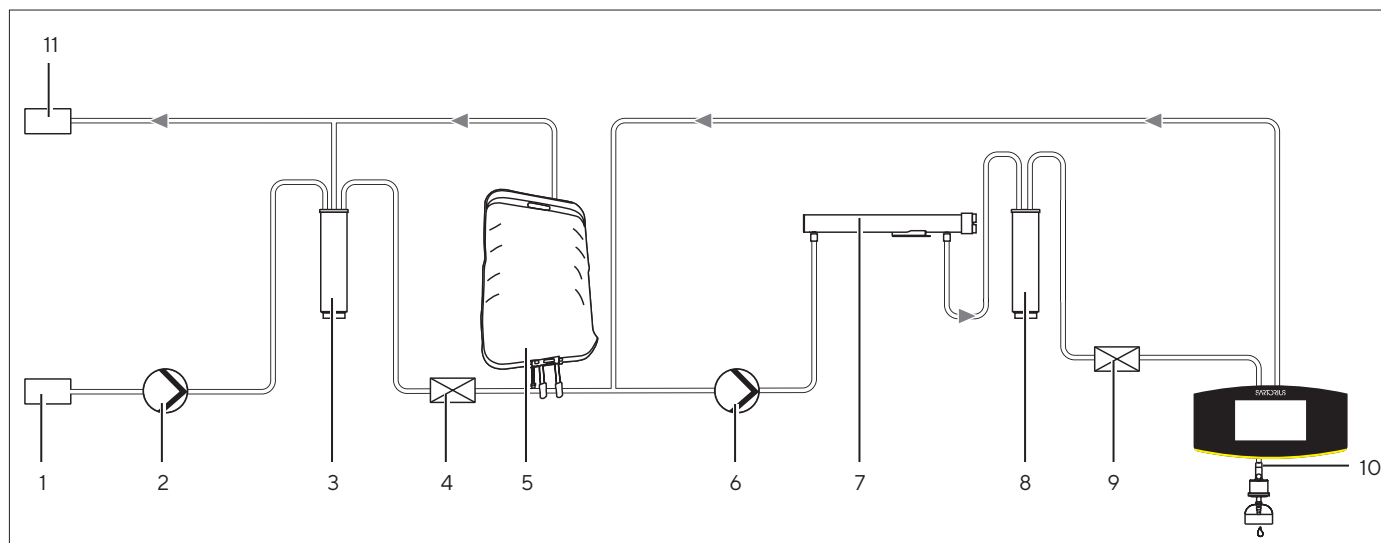


Fig.5: Arium® Mini Plus system setup

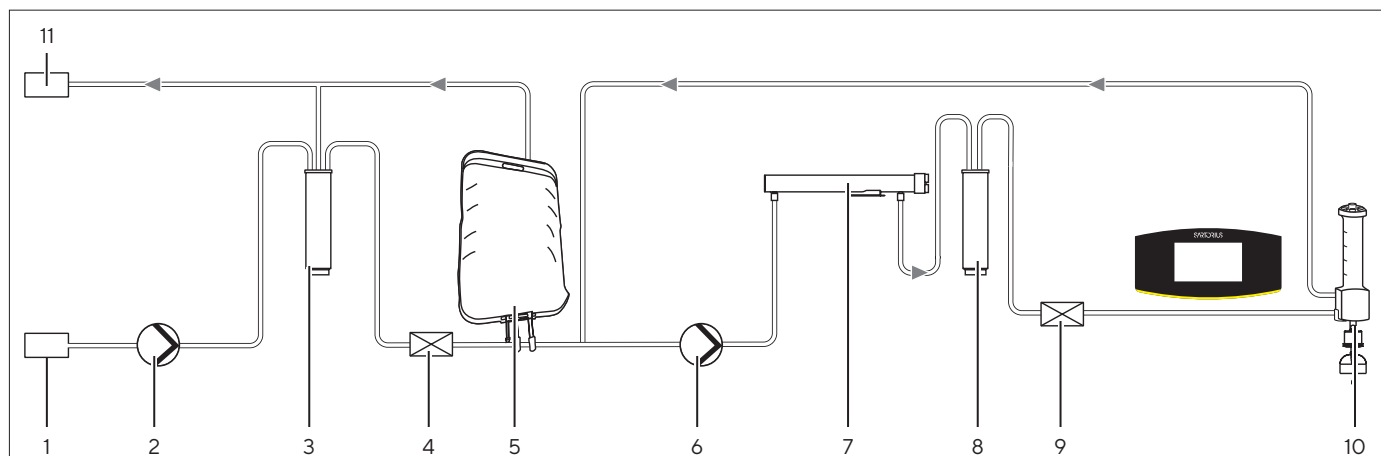


Fig.6: Arium® Mini Plus | Extend system setup

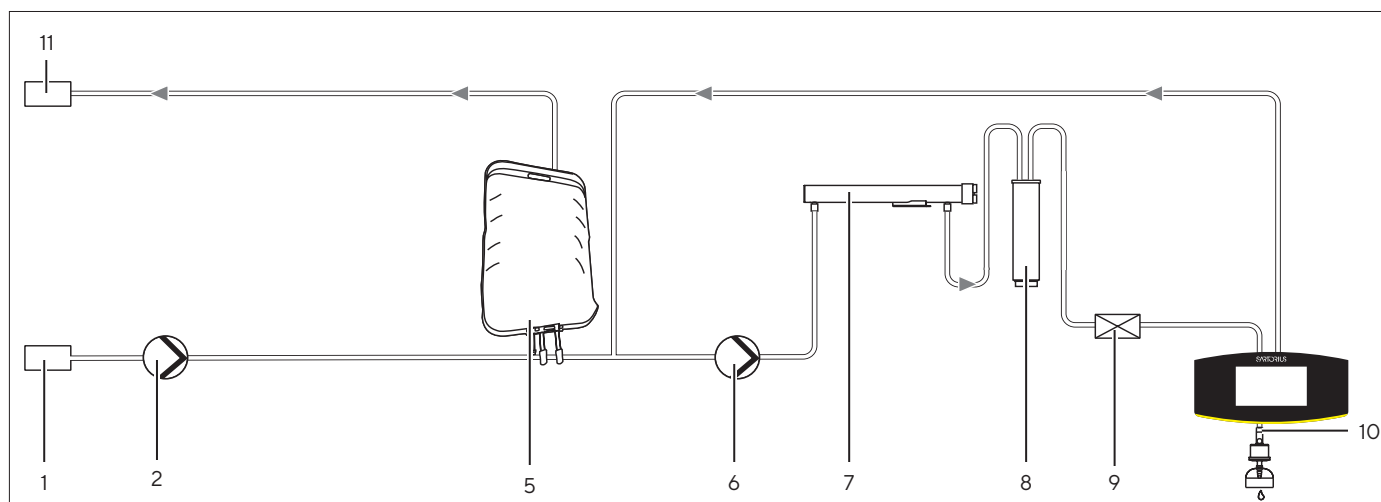


Fig.7: Arium® Mini system setup

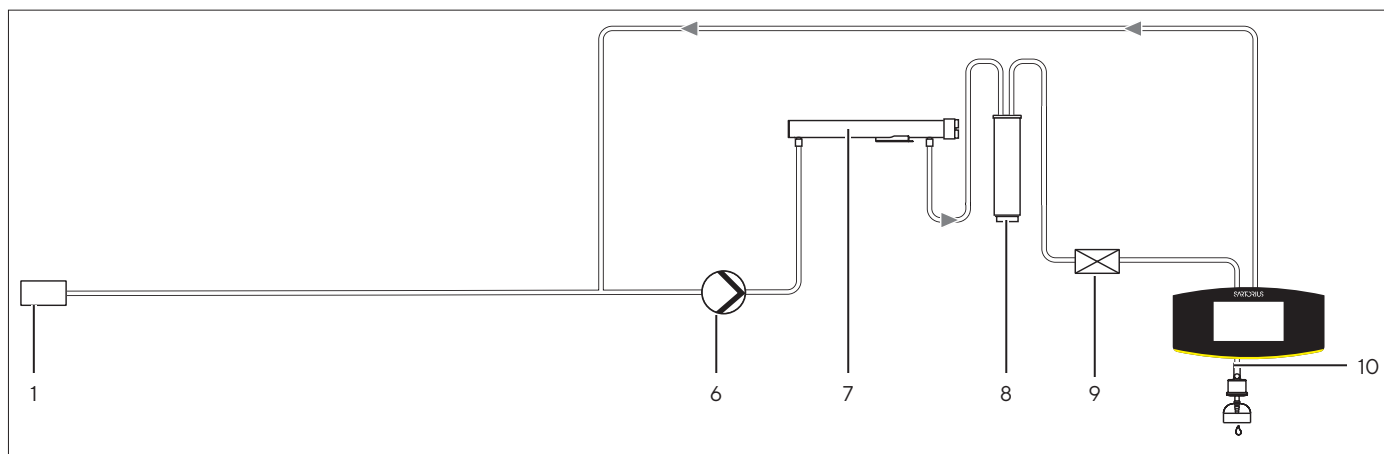


Fig. 8: Arium® Mini Essential system setup

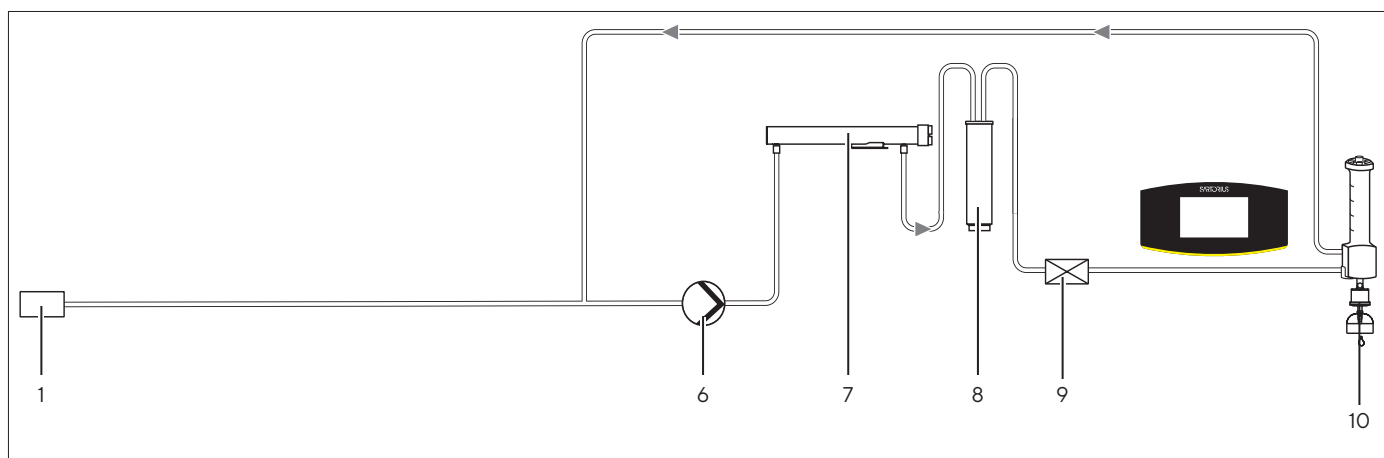


Fig. 9: Arium® Mini Essential | Extend system setup

No.	Description	No.	Description
1	Inlet (feed water)	7	UV lamp (185/254 nm)
2	Pump	8	Ultrapure water cartridge (Scientific Pack)
3	Pretreatment cartridge	9	Conductivity measurement
4	Conductivity measurement	10	Product water outlet
5	Bag	11	Outlet (drain water)
6	Pump		

3.5.2 Feed Water Intake

The feed water intake differs according to the device type:

Device type	Feed water intake
Arium® Mini Plus	Direct connection to tap water connection possible, automatic filling of the bag
Arium® Mini Plus Extend	Direct connection to tap water connection possible, automatic filling of the bag
Arium® Mini	Manual filling of the bag with pretreated water
Arium® Mini Essential	Direct connection to pretreated water
Arium® Mini Essential Extend	Direct connection to pretreated water

3.5.3 Treatment Stages

Device type	Required treatment stages
Arium® Mini Plus	Pretreatment stage, ultrapure water stage
Arium® Mini Plus Extend	Pretreatment stage, ultrapure water stage
Arium® Mini	Ultrapure water stage
Arium® Mini Essential	Ultrapure water stage
Arium® Mini Essential Extend	Ultrapure water stage

Pretreatment stage (first treatment stage)

In the first treatment stage, the tap water is treated with pretreated water: The feed water is passed through the pretreatment cartridge by means of a membrane pump. The pretreatment cartridge contains a combination of activated carbon, catalyst and downstream reverse osmosis. The pretreatment cartridge removes particles, salts and impurities, e.g. chlorine, from the tap water.

A large proportion of the impurities are discarded by means of a concentrate outlet (drainage water outlet).

The treated water is stored in the bag for further use. The quality of the treated water is monitored via a conductivity measuring cell.

Ultrapure water stage (second purification stage)

In the second purification stage, the pretreated water is purified to become ultrapure water (ASTM type 1): The pretreated water stored in the bag (Arium® Mini Plus with or without dispensing unit and Arium® Mini) or directly fed in pretreated water (Arium® Mini Essential with or without dispensing unit) is supplied to the ultrapure water circulation for this purpose.

The water fed in can be freed from organic residues using an optional UV lamp (185/254 nm). The pretreated water is then treated using the ultrapure water cartridge (Scientific Pack).

The quality of the ultrapure water is monitored via a conductivity measuring cell.

3.5.4 Final Purification Stage

The ultrapure water passes through a final filter as it is dispensed. A sterile filter or an ultrafilter can be used as a final filter.

3.5.5 Circulation of the Ultrapure Water

To guarantee ultrapure water of consistent and high quality, the ultrapure water purification system is constructed like a circulation ring:

- If **no** water is dispensed, the water circulates through the UV lamp and the ultrapure water cartridge.
- If there is **no** interaction with the device, after a short time the device switches automatically into ECO mode (see Chapter “7.6.1 Automatic ECO Mode,” page 58).

4 Operating Design

4.1 Menu

All system settings and work steps for the maintenance of the device can be carried out in the menu.

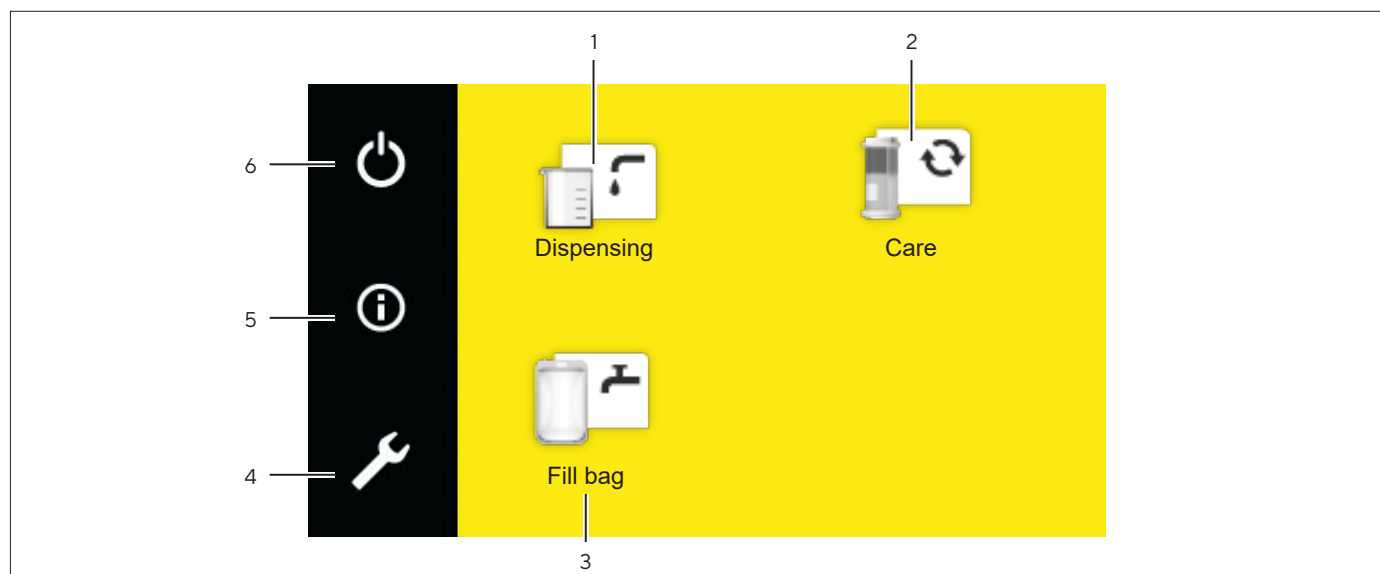








Fig.10: Menu in Arium® Mini (Example)

Pos.	Symbol	Name	Description
1		Dispensing	Opens the dispensing screen.
2		Care	Opens the "Care" menu.
3		Fill bag	Opens the wizard for bag filling (only with Arium® Mini).
4		Settings	Opens the "Settings" menu.
5		Information	Opens the "Information" menu.
6		Standby	Switches to standby mode.

4.2 Dispensing Mode

In dispensing mode, the display shows information about the water quality and the fill level of the bag, as well as the buttons relevant for dispensing. In the case of devices with a dispensing unit, removal can also be controlled via the hand-held part.

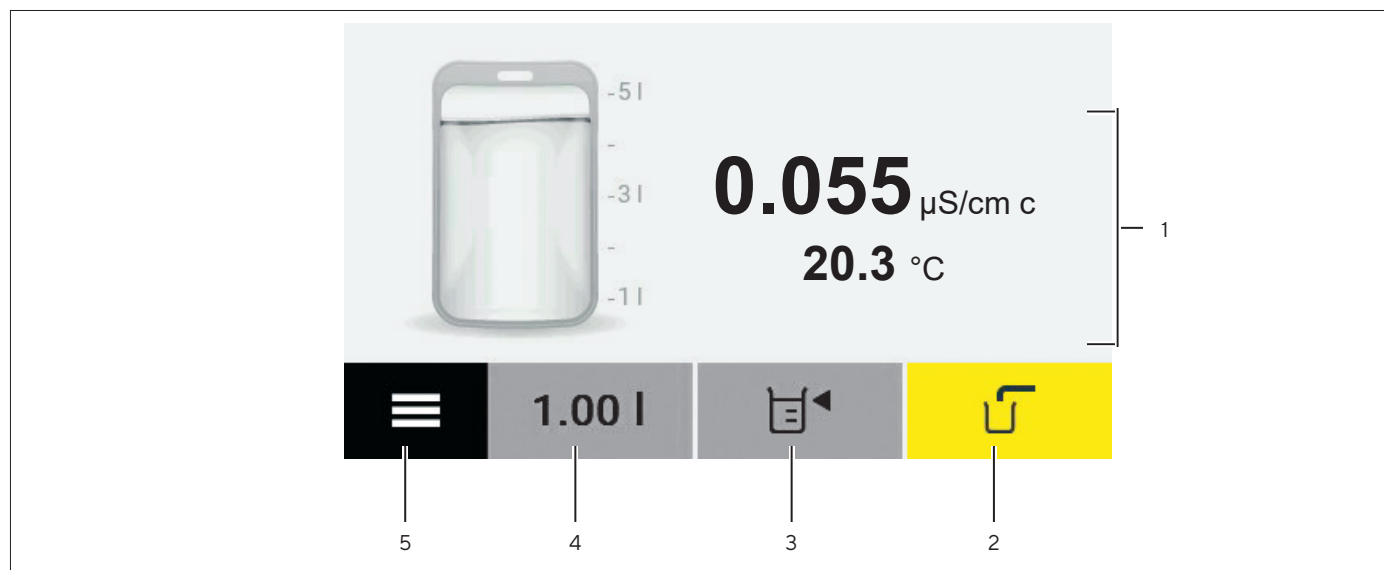


Fig.11: Dispensing mode Arium® Mini Plus (Example)

Pos.	Symbol	Name	Description
1		Operating area	Can display the following information: <ul style="list-style-type: none"> – Current ultrapure water conductivity – Current ultrapure water temperature – Fill level of the bag (only Arium® Mini Plus with or without dispensing unit and Arium® Mini) – Messages, warnings, errors
2		Manual dispense	Starts dispensing without preset dispense volumes.
3		Volume input	Opens a display for entering the dispense volume and starts the dispensing.
4		Volume-controlled dispensing	Starts dispensing with the displayed dispensing volume.
5		Menu	Opens the menu.

4.3 Message Display

The device displays three types of messages:

- Error messages (error):
 - Dispense is **not** possible.
 - The user only has limited options for troubleshooting errors.
- Warning messages (warning):
 - Dispense is possible.
 - The user can perform troubleshooting of warning messages.
- Status messages (info):
 - Dispense is possible.
 - There is information which the user should be aware of. Urgent action is **not** necessary.

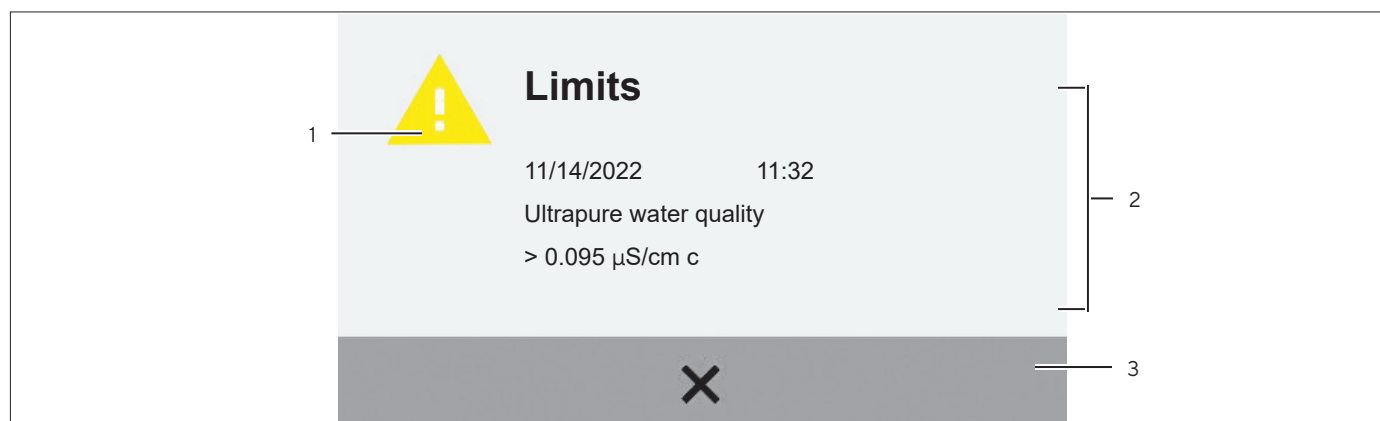






Fig.12: Warning message (Example)

Pos.	Symbol	Name	Description
1	  	Message symbol	Indicates which type of message is present: <ul style="list-style-type: none"> – Red symbol: Error message – Yellow symbol: Warning message – Gray symbol: Status message (info)
2		Message	Provides information about the message: <ul style="list-style-type: none"> – Error number or brief description of the message – Date and time of the occurrence – Message details
3		Acknowledge	Acknowledges the message.

4.4 Messages in Dispensing Mode

If several messages are active, the message list can be called up in the dispensing mode (see Chapter “4.5 Message List,” page 22). The message list only appears when several messages are active and at least one message **cannot** be displayed in the display for the conductivity or temperature of the water. If only one message is active, the message is opened directly instead of the message list.

If an error message is active, **no** dispensing is possible.

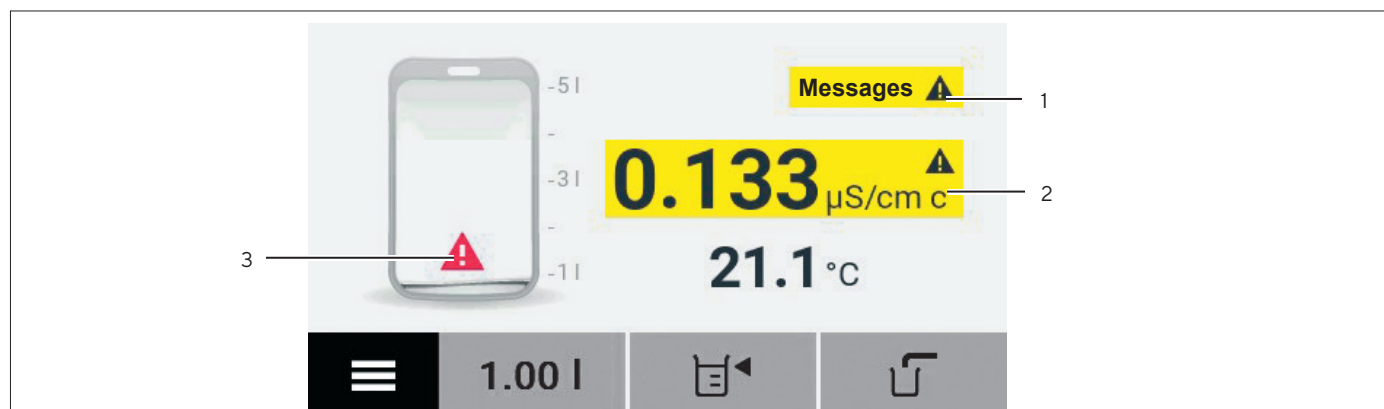


Fig.13: Message display in the dispensing mode (example)

Pos.	Symbol	Name	Description
1	 	Message list	Indicates that there is a message and opens the message list: <ul style="list-style-type: none"> – Red symbol: Error message – Yellow symbol: Warning message – Gray symbol: Status message
2	 	Value message	Indicates that there is a message about a value and opens the message list: <ul style="list-style-type: none"> – Red symbol: Error message – Yellow symbol: Warning message
3	 	Bag fill level message	Only for Arium® Mini Plus with or without dispensing unit and Arium® Mini: <ul style="list-style-type: none"> – Indicates the fill level of the bag – Red symbol: The bag is empty. Water dispense is not possible. – Yellow symbol: Only a little water can be dispensed.

4.5 Message List

All active messages can be viewed in the message list.

The messages in the message list are sorted according to priority. Error messages appear at the top. Within the same priority level, the messages are sorted according to date and time.

Messages **cannot** be deleted manually. They remain in the message list and will be shown on the display until their cause has been corrected. The device detects if the cause of a message has been corrected and automatically deletes the message from the message list and the display.

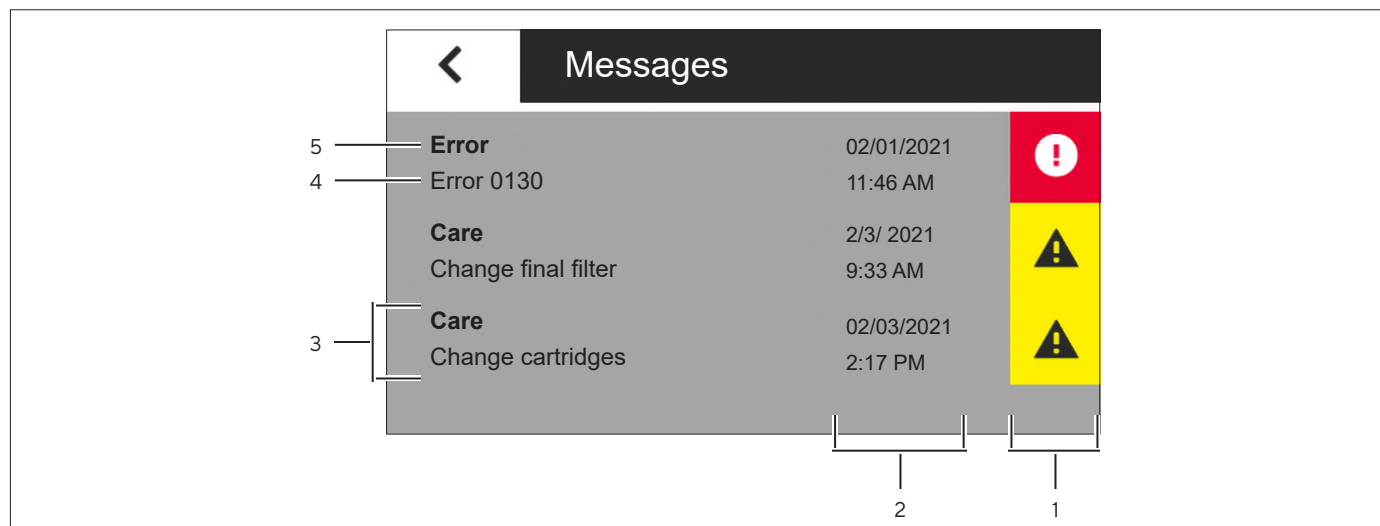





Fig.14: Message list (example)

Pos.	Symbol	Name	Description
1	  	Message symbol	Indicates which type of message is present: <ul style="list-style-type: none"> – Red symbol: Error Message – Yellow symbol: Warning Message – Gray symbol: Status message
2		Occurrence of the message	Displays the date and time when the message occurred.
3		Message	Displays and opens the message.
4		Brief description	Displays the error number or a short description of the message.
5		Message category	Displays the message category: <ul style="list-style-type: none"> – Error: Error message – Care: Message about consumables – Service: Message about device service – Quality: Message about the conductivity measurement – Temperature: Message about the temperature measurement – Info: Status message

4.6 Numeric Keypad

The numeric keypad is used to enter a dispensing volume or various system settings.

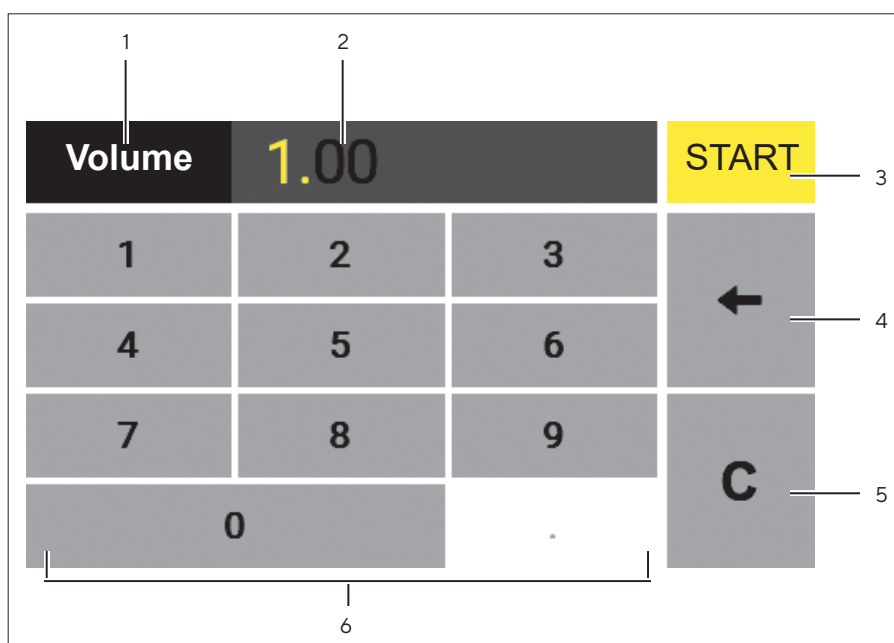





Fig.15: Volume input (example)

Pos.	Symbol	Name	Description
1		Name of the dialog box	Displays the name of the current dialog box.
2		Numerical value	Displays the currently entered numeric value.
3		Confirmation	Adopts the entered digit sequence and returns to the menu or to dispense mode.
4		Correction	Deletes the last digit entered.
5		Delete	Deletes all digits entered.
6		Numeric Keypad	Transfers numerical values to the dialog box.

Tip

When entering numeric values, it is only ever possible to select figures that are permitted for the value.

Example: Entering a dispensing volume of more than 5 liters is **not** possible. The number fields 6 to 9 are therefore inactive for this entry.

► Pay attention to the validity of the numeric values.

4.7 Volume Selection (Only for Devices With a Dispensing Unit)

The volume selection is intended to define the dispensing volume with the dispensing unit.

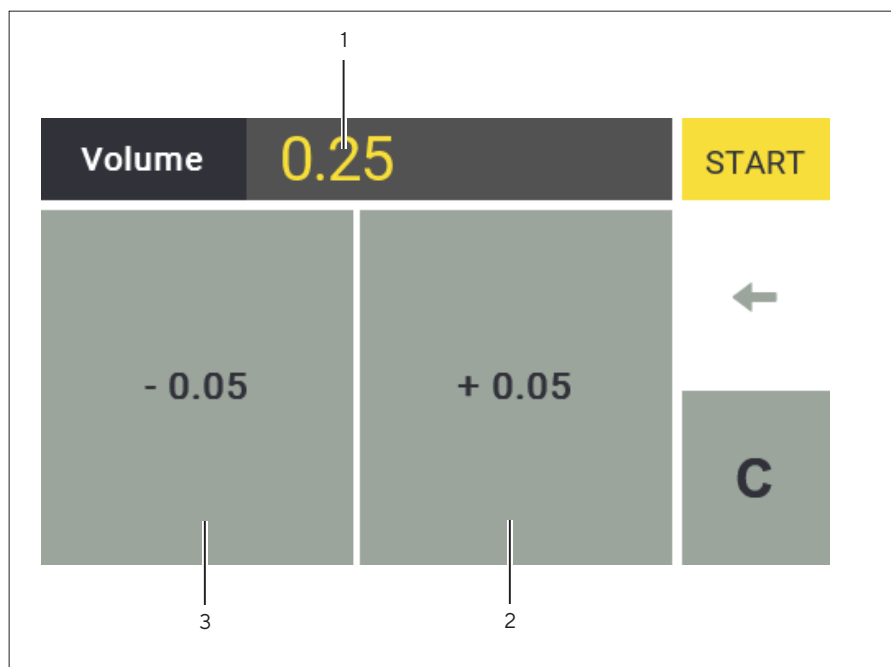


Fig.16: Volume for dispensing with the dispensing unit (example)

Pos.	Name	Description
1	Numerical value	Displays the currently entered numeric value.
2	Increase volume	Increases the numerical value by the displayed amount.
3	Decrease volume	Decreases the numerical value by the displayed amount.

4.8 Navigating the Menus

A display with touch function is provided for operating the device. If multiple menu items are available, you can use the touch screen to scroll up and down in order to select the desired entry.

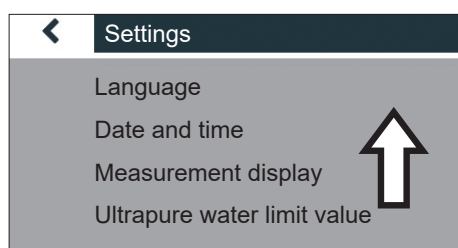
NOTICE

Sharp or pointed instruments such as ballpoint pens can damage the device!

Using unsuitable objects on the touch screen can damage it.

- ▶ The touch screen should only be operated by lightly pressing it using the tips of your fingers.
- ▶ The touch screen can also be operated while wearing laboratory gloves.

Procedure



- ▶ To scroll: Slowly swipe the touch screen up or down.
- ▷ The menu entries move in the corresponding direction.
- ▷ During scrolling, a gray scrollbar is displayed on the right for orientation.
- ▶ To select a menu item: Press the desired menu item.



- ▶ To confirm a selection: Press the [Confirmation] button.



- ▶ To cancel a process and return to the previous display: Press the [Cancel] button.



- ▶ To switch to the next highest menu level or to go back to the previous dialog box: Press the [Back] button.



- ▶ To go to manual value entry via the numeric keypad: Press the [Enter] button.



- ▶ To confirm a message: Press the [OK] button.






- ▶ To start a procedure: Press the [START] button.

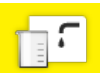




- ▶ To abort a process: Press the [STOP] button.

4.9 Menu Structure

► Navigating the menus (see Chapter “4.8 Navigating the Menus,” page 25).

Menu	Menu	Submenu	Description
 Standby			Starts standby mode.
 Information	Device Information		Shows all the features of the device, e.g. the model and the serial number.
	Measured values		Displays the current water quality from the ultrapure water stage and the pretreatment stage.
	Service information		If configured: Displays the Sartorius contact details and the next maintenance date.
	Reminder		Displays the next interval for replacing components, e.g., ultrapure water cartridge.
 Settings	Language		Changes the language of the operating interface.
	Date and time		Changes the date and time.
	Measurement display		Changes the units of the display for water quality and water temperature.
	Ultrapure water limit value		Specifies a limit value for ultrapure water quality.
	Final filter reminder		Activates deactivates the reminder for the next required final filter replacement. Allows for a selection of final filter types.
	Acoustic signals	Operating display button signals	Activates deactivates beeps when pressing a button.
		Key signals on hand-held part	Activates deactivates beeps when pressing the hand-held part keys [Plus], [Minus], and [Volume].
		Warning	Activates deactivates persistent beeps for warnings.
		Error	Activates deactivates beeps for error messages.
		Confirmation	Activates deactivates beeps for expired waiting times.
	Display brightness		Changes the brightness of the display.
	Volume adjustment		Readjusts the flow sensor of the device.
	Service mode		Only accessible to the Sartorius Service. Password-protected.
	Reset settings		Resets the device to default settings.

Menu	Menu	Submenu	Description
	Dispense		Opens the dispensing screen.
	Reminder		Shows when consumables next need to be replaced.
Care	Replace consumables	Bag	Starts a bag change (only Arium® Mini Plus with or without dispensing unit and Arium® Mini).
		Pretreatment cartridge	Starts a replacement of the pretreatment cartridge (only Arium® Mini Plus with or without dispensing unit).
		Ultrapure water cartridge	Starts a replacement of the ultrapure water cartridge.
		UV lamp	Starts a replacement of the UV lamp.
		Sterile filter / ultrafilter	Starts a replacement of the sterile filter or ultrafilter.
	Depressurization		Starts reducing the pressure in the device.
	Venting		Starts a manual rinsing process.
	Fill bag		Starts a manual filling of the bag (only with Arium® Mini).

4.10 Parameters of the “Settings” Menu

Parameters	Setting values	Explanation
Language	English, German, French, Italian, Spanish, Portuguese, Polish, Russian, Japanese, Chinese	
Date format	DD.MM.YYYY	Day.Month.Year
	MM/DD/YYYY	Month/Day/Year
	YYYY-MM-DD (ISO)*	Year-Month-Day, as per ISO standard
Time format	24 h*	24-hour mode
	12 h (a.m./p.m.)	12-hour mode
Displayed values (water quality)	$\mu\text{S}/\text{cm c}^*$	Displays the water quality on the dispense screen in $\mu\text{S}/\text{cm}$ and compensates to 25°C.
	$\mu\text{S}/\text{cm}$	Displays the water quality on the dispense screen in $\mu\text{S}/\text{cm}$, as well as the water temperature.
	$\text{M}\Omega\text{cm c}$	Displays the water quality on the dispense screen in $\text{M}\Omega\text{cm}$ and compensates to 25°C.
	$\text{M}\Omega\text{cm}$	Displays the water quality on the dispense screen in $\text{M}\Omega\text{cm}$, as well as the water temperature.
Displayed values (temperature)	°C	Displays the water temperature on the dispense screen in degrees Celsius.
	°F	Displays the water temperature on the dispense screen in degrees Fahrenheit.
	Off*	Deactivates the water temperature display (can only be selected if the water quality displayed is compensated to 25 °C).
Ultrapure water limit value		
Activation	On*	Activates the limit value for the water quality. If the limit value is exceeded during dispensing, a warning message appears. Dispense is possible.
	Off	Deactivates the limit value for the water quality.
Limit		Specifies the limit value for the water quality
Lock dispense	On	Activates the lock dispense. If the limit value is exceeded during dispensing, an error message is displayed. Dispense is not possible. This setting is recommended for particularly critical applications.
	Off*	Deactivates the lock dispense. If the limit value is exceeded: Dispensing is still possible.

Parameters	Setting values	Explanation
Final filter reminder	Active*	Activates deactivates the reminder for the final filter replacement (see Chapter 8.6, page 70).
	Inactive	
	Sterile filter*	Is used to select the final filter type.
	Ultrafilter	
	Reminder [weeks]	Defines the time interval for the reminder. Set to 4 weeks by default.
Acoustic signals		
Operating display button signals	On	Activates deactivates short beeps when pressing a button.
	Off*	
Key signals on hand-held part	On*	Activates deactivates beeps when pressing the hand-held part keys [Plus], [Minus], and [Volume].
	Off	
Warning	On*	Activates deactivates the persistent beep when a warning message appears, until the message is confirmed.
	Off	
Error	On*	Activates deactivates the persistent beep when an error message appears, until the message is confirmed. Does not deactivate the error signal if there is a leak.
	Off	
Confirmation	On*	Activates deactivates the longer beep at the end of time sequences, e.g. after a volume-controlled dispensing or after filling the bag.
	Off	
Display brightness	Bright*	Sets the display brightness to 100%.
	Dark	Sets the display brightness to 60%.
Volume adjustment		See Chapter 7.8, page 59.
Reset settings	Yes, reset	Resets all system settings to the default setting, e.g. the ultrapure water limit value and the displayed values. Does not reset the data for the reminders to replace consumables.
	No	Cancels the resetting of the settings.

* Default settings

5 Installation

5.1 Equipment Supplied

	Quantity		
	Arium® Mini	Arium® Mini Plus	Arium® Mini Essential
Device, bench-top system model	1	1	1
Inlet water tube: ¼" (outer diameter, length 2.40 m) with reducing connector ⅜" to ¼" (transition adapter) and tube (⅜" outer diameter, length 0.05 m)	–	1	1
Waste water tube: ⅜" outer diameter, length 2.40 m	1	1	–
Sampling tube: ¼" outer diameter, length 2.40 m	1	1	1
Tube adapter for inlet water with integrated sieve: ½" internal thread, ⅜" outer diameter	–	1	1
Tube adapter for inlet water with integrated sieve: G ¾" (internal thread, ⅜" outer diameter) with thread adapter G ¾" to ¼"	–	1	1
Bag outlet tube: ¼" outer diameter, length 1.50 m; with shut-off valve and tube (¼" outer diameter, length 0.10 m)	1	1	–
Tank filling tube: ¼" outer diameter, length 2.40 m	1	–	–
Tubing release tool	1	1	1
Power supply unit with country-specific power cable	1	1	1
Operating instructions	1	1	1
QA certificate	1	1	1
Only for devices with a dispensing unit			
Hand-held part with connection tube	1	1	1
Bracket with screw	1	1	1
Clip	1	1	1

Upon delivery, the following parts are located behind the left side cover of the device:

- Tubing
- Tubing adapter
- Hose release tool
- Power cord

- Only for devices with dispensing unit:
 - Hand-held part with connection tube
 - Bracket and screw
 - Clip

Consumables

The consumables are **not** included in the scope of delivery:

- Pretreatment cartridge
- Ultrapure water cartridge (Scientific Pack)
- Bag
- Final filter

5.2 Prerequisites for Installation at the Installation Location

Procedure

- Make sure that the following conditions are met at the place of installation:

Requirement	Features
Ambient conditions	<ul style="list-style-type: none"> – Suitability tested (ambient conditions, see Chapter 14.2, page 83; electromagnetic compatibility, see Chapter 14.4, page 84)
Footprint	<ul style="list-style-type: none"> – Stable and level surface – Sufficient space for the device (device space requirements, see Chapter “14.7 Device Properties,” page 88) – Sufficient load capacity for the device, including when it has been filled (device weight, see Chapter “14.7 Device Properties,” page 88)
Access to utilities	<ul style="list-style-type: none"> – Access to utilities is a maximum of 2 meters: <ul style="list-style-type: none"> – Feed water supply – Power outlet – Pressureless drain
Feed water quality	<ul style="list-style-type: none"> – Suitability tested (see Chapter “14.6 Feed Water Quality,” page 86)

5.3 Unpacking and Setting up the Device

We recommend that installation of the device be carried out by Sartorius Service. Contact Sartorius Service in this regard.

CAUTION

Risk of electric shock due to leaking water!

Water may spill when using the device. Electric shocks can occur if water comes into contact with electrical appliances.

- ▶ Do **not** place the device close to electrically powered devices.


CAUTION

Danger of fire or explosion!

The device contains components that can ignite highly inflammable or combustible materials.

- ▶ Do **not** operate the device in the vicinity of such substances.

Procedure

- ▶ If the device is stored temporarily: Observe the storage information (see Chapter "10.1 Storage," page 79).
- ▶  **CAUTION** Injury to the feet due to falling device components during improper transport!
 - ▶ Do **not** lift or transport the device by the side covers or components of the dispensing unit.
 - ▶ Hold the device underneath the display on the front side and by the recess for the power supply unit on the back side and lift it carefully.
- ▶ Take the device out of the packaging and set it up at the intended installation site.



5.3.1 Removing the Left Side Cover

Procedure

- ▶ Reach behind the side cover and pull it to one side (1).
- ▶ Pull the side cover upwards and remove (2).



5.3.2 Removing Parts

Procedure

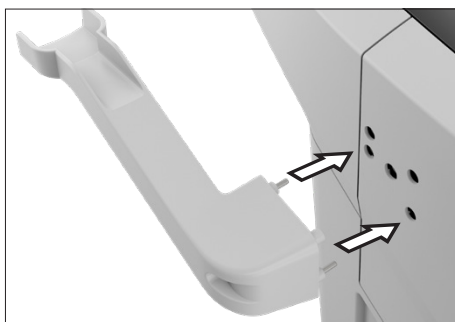
- ▶ Remove the following parts from inside the device:
 - Tubing
 - Tube adapter
 - Tubing release tool
 - Power cable
- ▶ Only for devices with dispensing unit: Remove the following parts from inside the device:
 - Hand-held part with connection tube
 - Bracket and screw
 - Clip

5.4 Mounting the Dispensing Unit

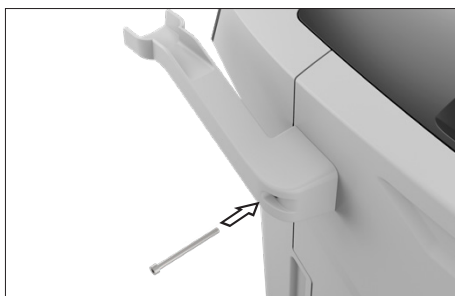
Tool

Allen key, 1 piece

Procedure



- ▶ Place the protruding pins of the brackets in the recesses provided on the side of the device.



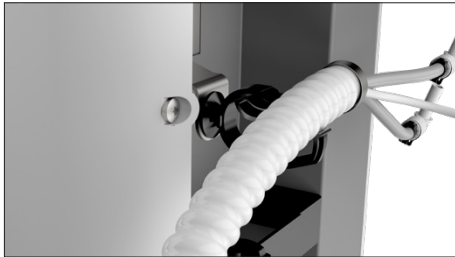
- ▶ Secure the bracket to the device using the screw.



- ▶ Place the hand-held part in the magnetic bracket.



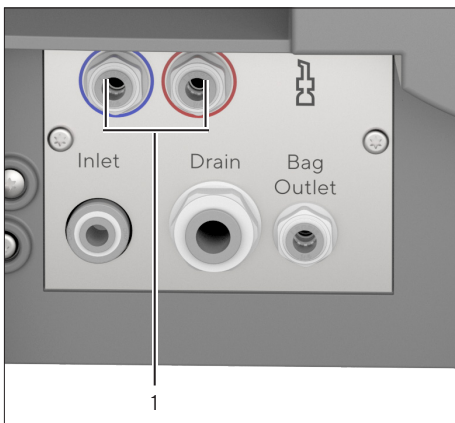
- ▶ Press the clip into the plate on the back of the device.
- ▷ The clip engages.
- ▶ Open the clip.



- ▶ **⚠ CAUTION** Grazing to the hand due to the edges of the protruding plate. Avoid the plate when attaching the connection tube.
- ▶ Press the connection tube into the clip.
- ▷ The clip closes around the connection tube.
- ▶ Press to secure the clip.

5.4.1 Connecting the Tube Unit

Procedure



- ▶ Connect the color-marked feed water tube (blue) and return water tube (red) of the dispensing unit to the identically marked connections (1).
- ▶ **⚠ CAUTION** Risk of injury due to liquid spurting out. Check the secure position of the connection tube.

6 Startup

We recommend having the initial startup of the device carried out by Sartorius Service. Please contact Sartorius Service in this regard.

The startup steps required are shown on the device display.

6.1 Connecting the AC Adapter

CAUTION

Risk of electric shock due to incorrect handling of power cables!

The use of damaged or **non-standard** power cables as well as the mishandling of power cords can cause fatal electric shock or equipment damage.

- ▶ Only connect the device (protection class 1) to properly installed power sockets with protective grounding conductors (PE) with a fuse of a maximum of 16 A.
- ▶ Do **not** replace the removable power cable with a power cable with insufficient nominal value or of the wrong type.
- ▶ **Never** disconnect the device from the protective grounding conductor.
- ▶ Connect to the power supply according to the regulations of your specific country.
- ▶ **Never** plug the power cable into the mains wall outlet when it is disconnected from the device.
- ▶ Make sure that the power plug or another suitable disconnecting device for the power can be easily reached in case of danger.

NOTICE

Equipment damage due to operation with third-party equipment!

The use of third-party power supplies **not** authorized by Sartorius may cause damage to the device.

- ▶ Use only the original Sartorius power supply.

Procedure

- ▶ Check whether the plug design of the power cord is equivalent to your country's standard.
 - ▶ If required: Call Sartorius Service or your dealer.
- ▶ Connect the power supply to the connection labeled "Power" on the rear of the device.





- ▶ Fix the power supply to the housing using the supplied Velcro strip. To do this, pull the Velcro strip through the eyelet as shown in the figure.
- ▶ Lay the protruding power cord so that it **cannot** be damaged or obstruct subsequent connection of hoses.

Connecting the Dispensing Unit

Procedure



- ▶ Connect the connection cable of the hand-held part to the dispensing unit connection (1).
- ▶ Use the power cable to connect the device to the power supply.
- ▶ The device starts and performs a system check.

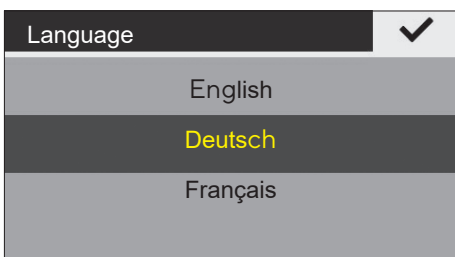
6.2 Configuring Device Settings (Startup)

6.2.1 Setting the Language

Once the system check is complete, the “Language” dialog box appears. The wizard automatically carries out all of the startup steps. Startup of the device takes up to 120 minutes and **cannot** be interrupted.

All the system settings (e.g. date, time, displayed values) configured during startup can be changed later in the “Settings” menu (see Chapter “7.7 Changing System Settings,” page 58).

Procedure



- ▶ Select the language for the display text. The factory setting is English.
- ▶ Confirm the selection with the [Confirmation] button.
- ▶ The “Startup” dialog box is displayed.

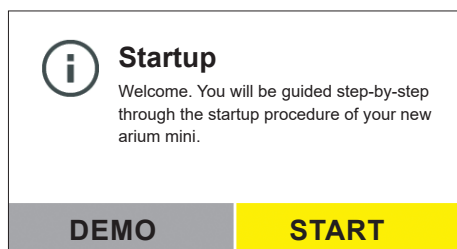
6.2.2 Launching Startup Mode

Requirements

The “Startup” dialog box is displayed.

Procedure

- ▶ To launch startup mode: Press the [START] button.
- ▶ To launch demo mode: Press the [DEMO] button. Demo mode is only accessible by Sartorius employees and is password-protected.



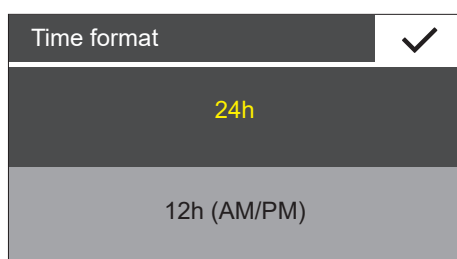
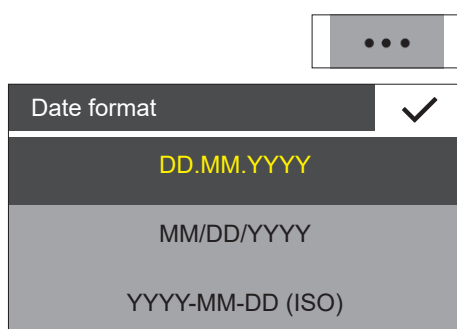
6.2.3 Setting the Date and Time

Requirements

The “Date and time” dialog box is displayed.

Procedure

- ▶ Press the [Input] button.
- ▶ Select the desired date format (for setting options see Chapter “4.10 Parameters of the “Settings” Menu,” page 28).
- ▶ Confirm the selection with the [Confirmation] button.
- ▶ Type the date into the entry field.
- ▶ Confirm the input with the [Confirmation] button.



- ▶ Select the desired time format (for Setting options, see Chapter “4.10 Parameters of the “Settings” Menu,” page 28).
- ▶ Confirm the selection.
- ▶ Type the time into the entry field.
- ▶ Confirm the input with the [Confirmation] button.

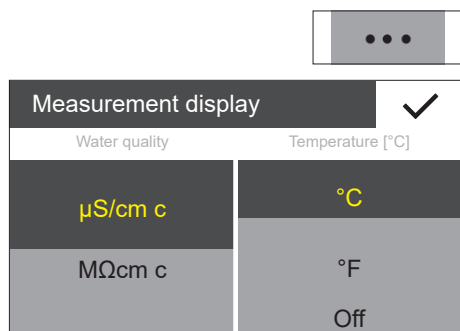
6.2.4 Setting the Displayed Values

Requirements

The “Displayed values” dialog box appears.

Procedure

- ▶ Press the [Input] button.
- ▶ Set the desired format for the water quality (for Settings options, see Chapter “4.10 Parameters of the “Settings” Menu,” page 28).
- ▶ Confirm the selection with the [Confirmation] button.
- ▶ Set the desired format for the temperature (for Settings options, see Chapter “4.10 Parameters of the “Settings” Menu,” page 28).
- ▶ Confirm the input with the [Confirmation] button.



6.2.5 Concluding Device Settings (Startup)

Requirements

The “Remove cover” dialog box is displayed. Depending on the type of device, an ultrapure water cartridge and | or pretreatment cartridge needs to be inserted.

Procedure

- ▶ Insert the required cartridge (see Chapter “6.3 Inserting Pretreatment Cartridge (only Arium® Mini Plus),” page 39 or Chapter “6.4 Inserting Ultrapure Water Cartridge,” page 40).

6.3 Inserting Pretreatment Cartridge (only Arium® Mini Plus)

A pretreatment cartridge must be inserted in the device. The cartridge holder for the pretreatment cartridge is marked "R".

Procedure

- ▶ Remove the front cover.
- ▶ Confirm removal of the cover with the [OK] button.
- ▷ The dialog box "Insert cartridge" appears.
- ▶ **NOTICE** Equipment damage due to impurities! If dirt or foreign objects enter the ultrapure water circulation, individual device components can become clogged or wear out faster. Do not **touch** the connections of the pretreatment cartridge. This will prevent impurities from getting into the device.
- ▶ Remove the plugs from all three connections of the pretreatment cartridge.
- ▶ Hold the pretreatment cartridge in the upper area and insert it into the holder marked "R". To do this, push the pretreatment cartridge straight into the guide rails of the cartridge holder.



- ▶ Press the pretreatment cartridge firmly into the cartridge holder until it locks into the guide rails with a distinct clicking sound on both the left and right sides. Check that the pretreatment cartridge has been installed securely by turning it slightly.
- ▶ Confirm the insertion of the pretreatment cartridge with the [OK] button.
- ▷ The "Insert cartridge" dialog box (Step 2) appears. An ultrapure water cartridge must be inserted (see Chapter 6.4, page 40).

6.4 Inserting Ultrapure Water Cartridge

An ultrapure water cartridge must be inserted in the device. The cartridge holder for the ultrapure water cartridge is marked "L".

Procedure

- ▶ Remove the front cover.
- ▶ Confirm removal of the cover with the [OK] button.
- ▷ The dialog box "Insert cartridge" appears.
- ▶ **NOTICE** Equipment damage due to impurities! If dirt or foreign objects enter the ultrapure water circulation, individual device components can become clogged or wear out faster. Do **not** touch the connections of the ultrapure water cartridge. This will prevent impurities from getting into the device.
- ▶ Remove the plugs from the two external ultrapure water cartridge connections.
- ▶ Firmly hold the upper area of the ultrapure water cartridge and insert into the holder marked "L". In order to do this, push the ultrapure water cartridge straight into the guide rails of the cartridge bracket.



- ▶ Press the ultrapure water cartridge firmly into the cartridge holder until it locks into the guide rails with a distinct clicking sound on both the left and right sides. Check that the ultrapure water cartridge has been installed securely by turning it slightly.
- ▶ Confirm the insertion of the ultrapure water cartridge with the [OK] button.

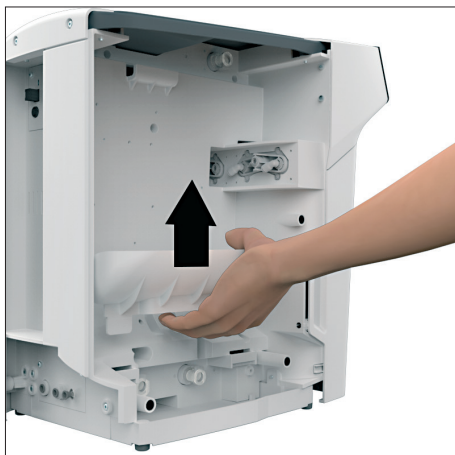


- ▷ The dialog box "Attaching the cover" appears.
- ▶ Put the front cover on the device.
- ▶ Confirm that the cover has been put on with the [OK] button.
- ▷ If an Arium® Mini Plus with or without dispensing unit or Arium® Mini is used: The dialog box "Insert bag" appears. A bag must be inserted (see Chapter 6.5, page 41).
- ▷ If an Arium® Mini Plus with or without dispensing unit is used: The dialog box "Connect tubing" appears. Feed water tubing must be connected (see Chapter 6.6.1, page 42).

6.5 Inserting Bag (only Arium® Mini Plus or Arium® Mini)

Procedure

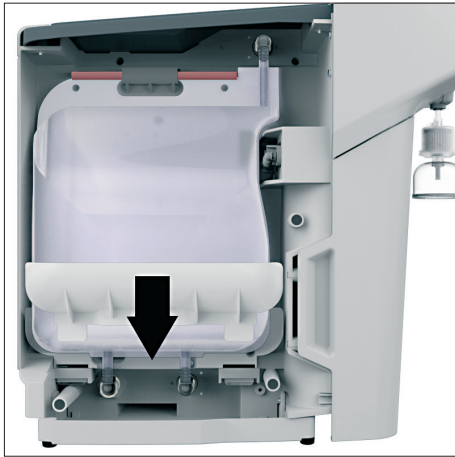
- ▶ Remove the left side cover (see Chapter “5.3.1 Removing the Left Side Cover,” page 32).
- ▶ To make it easier to reach the lower connections on the device: Pull the tray up and remove it from the device.



- ▶ **NOTICE** Danger of damage to the device from improper handling! The bag may be damaged if forced into the guide rails: The bag can be damaged. Use only slight pressure when inserting the bag into the guide rail.
- ▶ Using the handle, slide the bag into the guide rail in the upper portion of the device housing. The bag is secure when the reinforcing piece on the carrying handle is completely engaged in the guide rail.



- ▶ Using the quick connector, fasten the upper hose of the bag to the device's upper connection.
- ▷ The quick connector locks into place with an audible click.
- ▶ Fasten the two lower bag hoses to the lower device connections using the quick connectors.
- ▷ The quick connectors lock into place with an audible click.



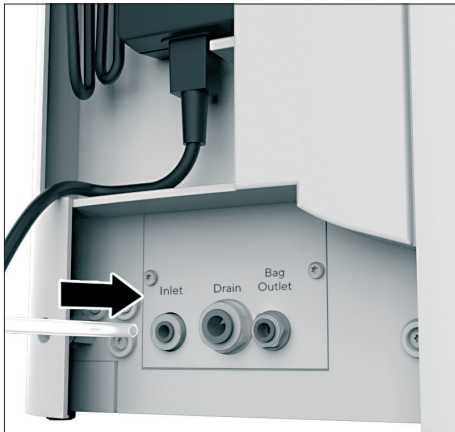
- ▶ Reinsert the tray. Make sure **not** to damage the bag.
- ▶ Reattach the left side cover of the device.
- ▶ Confirm the insertion of the bag with the [OK] button.
- ▶ The dialog box “Connect tubing” appears. Tubing must be connected (see Chapter 6.6, page 42).

6.6 Connecting Tubing

6.6.1 Connecting Feed Water Tubing (only Arium® Mini Plus or Arium® Mini)

To prevent the entry of particles into the device, it is recommended to use the supplied screens.

Procedure



- ▶ Connect the feed water tubing to the connection labeled “Inlet” on the rear of the device.
- ▶ **NOTICE** Water leakage due to excessive inlet pressure! If the water pressure is too high: Water can leak out of the device. Check the water pressure inlet specification (see Chapter “14.6 Feed Water Quality,” page 86).
 - ▶ If required: Reduce the inlet pressure.
- ▶ **NOTICE** Water leakage can occur if feed water tubing is not watertight! If the feed water tubing is deformed or **not** inserted deep enough: Water can leak out. After startup, make sure that all external water connections are leak proof.
- ▶ Connect the feed water tubing to the feed water supply.

6.6.2 Connecting Tank Filling Tubing (only Arium® Mini)

To prevent the entry of particles into the device, it is recommended to use the supplied screens.

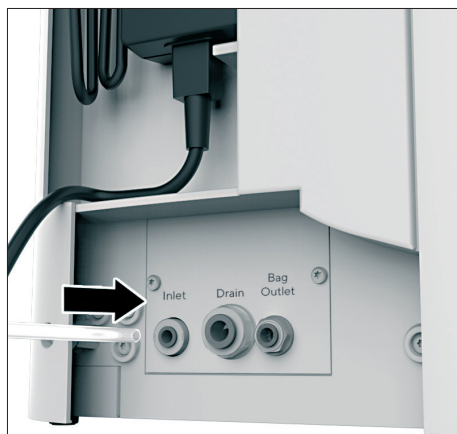
NOTICE

Contamination of tank filling tubing!

If the tank filling tubing is **improperly** stored: The tank filling tubing can be contaminated. Sufficient quality of the pure water that the bag is filled with is **no longer** guaranteed in this case.

- ▶ Do **not** leave the tank filling tubing on the floor or near sources of contamination (e.g. in the vicinity of an outlet).
- ▶ If the tank filling tubing **cannot** be kept secure while connected to the device: Remove the tank filling tubing and keep it in a clean place.
- ▶ If the tank filling tubing is contaminated:
 - Replace the bag (see Chapter 8.5.2, page 63).
 - Replace the tank filling tubing.
- ▶ Use the tank filling tubing exclusively for filling.

Procedure



- ▶ Connect the tank filling tubing to the connection labeled “Inlet” on the rear of the device.
- ▶ **NOTICE** Water leakage due to excessive inlet pressure! If the water pressure is too high: Water can leak out of the device. Check the water pressure inlet specification (see Chapter “14.6 Feed Water Quality,” page 86).
- ▶ **NOTICE** Water leakage can occur if feed water tubing is not watertight! If the feed water tubing is deformed or **not** inserted deep enough: Water can leak out. After startup, make sure that all external water connections are leak proof.

6.6.3 Connecting Drain Water Tubing (only Arium® Mini Plus or Arium® Mini)

NOTICE

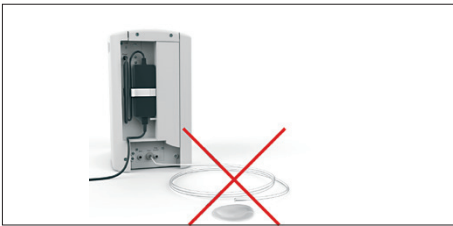
Bag can burst due to excess pressure!

If the drain water tubing of the device is sealed off, clogged or exposed to counter-pressure, the bag may burst.

- ▶ Do **not** seal off, clog or expose the drain water tubing to counter-pressure.

Procedure

- ▶ Connect the drain water tubing to the connector labeled “Drain” on the rear of the device.



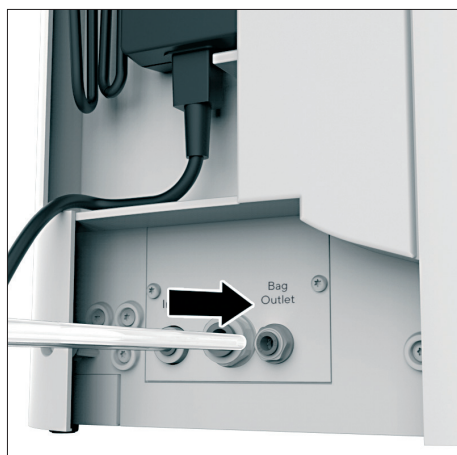
- ▶ **NOTICE** Water can leak out if drain water hose detaches! Water can escape from the drain water tubing during operation. Attach the drain water tubing to the outlet.



- ▶ Guide the free end of the waste water tubing to the drain and attach it.

6.6.4 Connecting the Tank Outlet Tubing (only Arium® Mini Plus or Arium® Mini)

Procedure



- ▶ Connect the long tank outlet tubing to the connector labeled “Bag Outlet” on the rear of the device.
- ▶ Close the ball cock.
- ▷ Later on in the process, the wizard will display a prompt for opening and closing the ball valve.

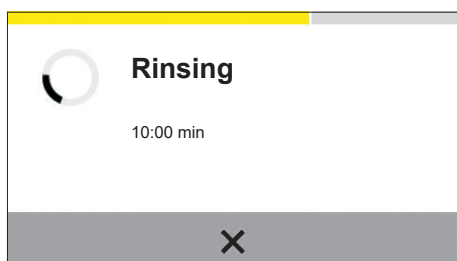
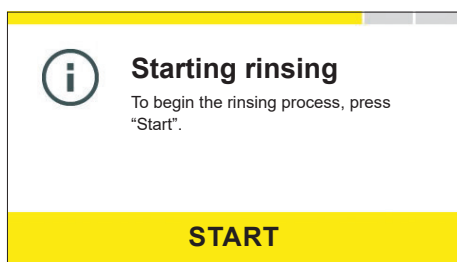
6.7 Rinsing Pretreatment Cartridge (only Arium® Mini Plus)

Requirements

The “Prepare rinsing” dialog box is displayed.

Procedure

- ▶ Open the ball valve on the bag outlet tube.
- ▶ Confirm the opening of the ball valve with the [OK] button.
- ▶ Guide the bag outlet tube towards the drain for rinsing. Water may leak out of the tube during rinsing.
- ▷ The “Start rinsing” dialog box is displayed.
- ▶ Start the rinsing process of the pretreatment cartridge. To do this, press the [START] button.



- ▷ During the rinsing process, the remaining rinse time is displayed in minutes.
- ▶ To interrupt the rinsing process:
 - ▶ Press the [Cancel] button.
 - ▷ The “Start rinsing” dialog box is displayed again.
- ▶ To resume the rinsing process after an interruption: Press the [START] button.
- ▶ When the rinsing process is complete:
 - ▷ The “Close bag outlet” dialog box is displayed.
 - ▶ Close the ball valve on the bag outlet tube.
 - ▶ Confirm the closing of the bag outlet. To do this, press the [OK] button.

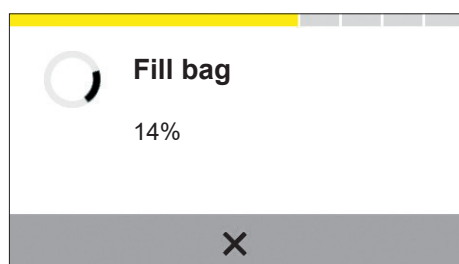
6.8 Rinsing the Ultrapure Water Cartridge

6.8.1 Filling of the Bag (only Arium® Mini Plus and Arium® Mini)

Automatic Filling (Arium® Mini Plus)

During rinsing, the device fills and rinses the ultrapure water cartridge. Air is removed from the ultrapure water circulation in the process.

- ▶ If the bag **no longer** contains a sufficient amount of water (Tank Fill Level) or empties during the rinsing process, the bag is filled automatically.
- ▶ If the fill level is sufficient to continue rinsing: The “Prepare rinsing” dialog box is displayed. It is necessary to connect a sampling tube (see Chapter “6.6.2 Connecting Tank Filling Tubing (only Arium® Mini),” page 43).



Manually fill the bag (only for Arium® Mini)

NOTICE

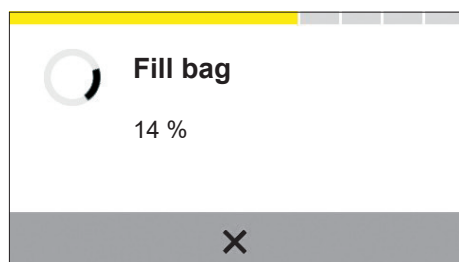
Air ingress may interfere with device functions!

If there is **not** enough water available during the filling process, air may be drawn in. Air in the system can interfere with the functioning of the device.

- ▶ Prepare a vessel containing a sufficient quantity of pretreated water (at least 5 liters).

Procedure

- ▶ Place the tank filling tube into the container and secure it to ensure that it is fully submerged throughout the filling process.
- ▶ Press the [START] button.
- ▶ The bag is filled with pretreated water. The progress is displayed as a percentage.
- ▶ To cancel the filling process early:
 - ▶ Press the [Cancel] button.
 - ▶ The “Fill bag” dialog box appears.
- ▶ To resume the filling process after an interruption: Press the [START] button.
- ▶ When the filling process is complete, the “Prepare rinsing” dialog box appears. It is necessary to connect a sampling tube (see Chapter “6.8.2 Performing a Rinse Cycle,” page 47).



6.8.2 Performing a Rinse Cycle

During rinsing, the device fills and rinses the ultrapure water cartridge. Air is removed from the ultrapure water circulation in the process.

Requirements

The “Prepare rinsing” dialog box is displayed.

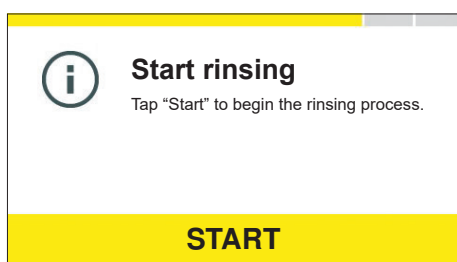
Procedure



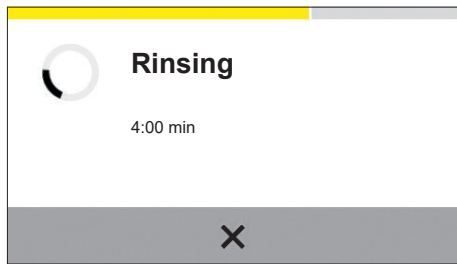
- For devices without a dispensing unit: Press the sampling tube into the quick connector of the water outlet.



- For devices with a dispensing unit: Press the sampling tube into the quick connector of the water outlet on the hand-held part.
- Guide the free end of the sampling tube to the drain or place it in a container (at least 6 liters).
- To confirm the connection of the sampling tube: Press the [OK] button.



- Start the rinse cycle for the ultrapure water cartridge. To do this, tap the [START] button.

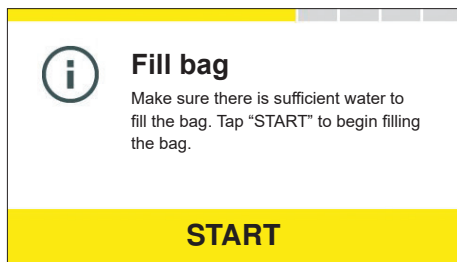


- ▷ During the rinse cycle, the remaining rinse time is displayed in minutes.
- ▶ To interrupt the rinse cycle:
 - ▶ Press the [Cancel] button.
 - ▷ The dialog box "Start rinsing" appears again.
- ▶ To resume the rinse cycle after an interruption: Tap the [START] button.
- ▷ The rinse cycle is performed.
- ▷ Once the rinsing process is complete: The "End rinsing" dialog box is displayed.
- ▶ Remove the sampling tubing.
- ▶ To confirm the removal of the sampling tube: Press the [OK] button.
- ▶ Connect the final filter (see Chapter "6.9 Connecting the Final Filter," page 49).

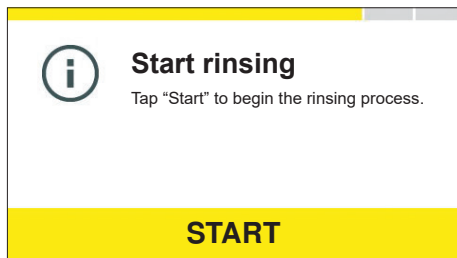
6.8.3 Re-filling the Bag During the Rinse Cycle (only Arium® Mini)

If the bag does **not** contain enough water or empties during the rinse cycle:
The bag must be refilled.

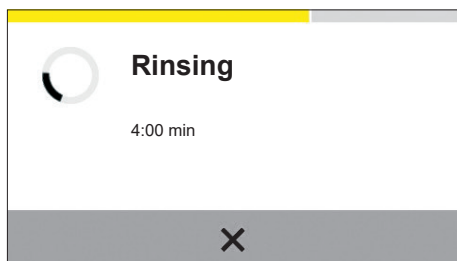
Procedure



- ▶ The dialog box "Fill bag" appears: Fill the bag (see Chapter 6.8, page 46).



- ▶ Start the rinse cycle for the ultrapure water cartridge. To do this, tap the [START] button.



- ▷ During the rinse cycle, the remaining rinse time is displayed in minutes.
- ▶ To interrupt the rinse cycle:
 - ▶ Press the [Cancel] button.
 - ▷ The dialog box "Start rinsing" appears again.
- ▶ To resume the rinse cycle after an interruption: Tap the [START] button.
- ▷ The rinse cycle is performed.

6.8.4 Completing the Rinsing Process

Procedure

- ▶ Once the rinsing process is complete and the “End rinsing” dialog box is displayed:
 - ▶ Remove the sampling tubing. To do this, press the dark ring of the quick connector upwards using the tubing release tool and pull out the sampling tube downwards. For devices with a dispensing unit, the water outlet is on the hand-held part.
 - ▶ Confirm the rinsing process via the [OK] button.
- ▷ The “Connect filter” dialog box appears. A final filter needs to be connected (see Chapter 6.9, page 49).



6.9 Connecting the Final Filter

Requirements

The “Connect filter” dialog box appears on the display.

Procedure

- ▶ Attach the bell assembly to the final filter.
- ▶ Press the final filter into the quick connector of the water outlet. For devices with a dispensing unit, the water outlet is on the hand-held part.
- ▶ Confirm that the final filter has been connected with the [OK] button.
- ▷ The display changes to the dispensing screen.



6.10 Rinsing the Final Filter

Requirements

The device is now in dispensing mode.

Procedure

- ▶ Place a vessel under the final filter.
- ▶ If a sterile filter is used as the final filter: Open the final filter's vent valve.
- ▶ Remove the protective cap from the bell assembly.
- ▶ If a sterile filter is used as the final filter: Remove at least 5 liters of water. This will rinse the final filter.
- ▶ If an ultrafilter is used as a final filter: Remove at least 20 liters of water. This will rinse the final filter.
- ▶ Only Arium® Mini Plus with or without dispensing unit or Arium® Mini: If the bag **no longer** contains sufficient water or empties during the rinsing process:
 - ▶ Only Arium® Mini: To fill the bag: Press the [YES] button. Fill the bag manually (see Chapter "7.5.2 Starting Manual Filling in Menu," page 56).
 - ▶ Only Arium® Mini Plus with or without dispensing unit: Wait until the bag has reached a sufficient filling level.
 - ▶ Finish the rinsing process.
- ▶ When the rinsing process is complete: Close the vent valve.
- ▶ Attach the protective cap to the bell assembly.
- ▷ Startup is complete.

7 Operation

7.1 Switching the Device On and Off

Procedure

- ▶ To switch the device on: Connect the device to the power supply.
- ▷ The device starts and performs a system check.
- ▶ To switch the device off: Disconnect the device from the power supply.

Tip

If the device is switched off in normal operation, e.g., in the evening or at weekends, then consistent ultrapure water quality is no longer guaranteed. To ensure a consistent ultrapure water quality, we recommend activating the standby mode (see Chapter “7.6 Activating or Deactivating Standby Mode,” page 58).

7.2 Dispensing Ultrapure Water

7.2.1 Preparing to Dispense Ultrapure Water

The ultrapure water can be dispensed as follows:

- Manual dispensing
- Volume input

Requirements

- The device is ready for operation (see “6 Startup”, page 35).
- The device is in dispensing mode.

NOTICE

Water leakage due to vessel overflow!

- ▶ Do not allow water to be dispensed unattended. The vessel being filled may overflow.
 - ▶ To avoid dispensing volumes that are too large: Dispense the ultrapure water in a volume-controlled manner (see Chapter 7.2.3, page 53).
-

Procedure

- ▶ If dispensing takes place via the final filter: Secure the protective cap on the filling bell of the final filter.
- ▶ Place a suitable vessel under the water outlet.
- ▶ The device is ready to dispense ultrapure water.

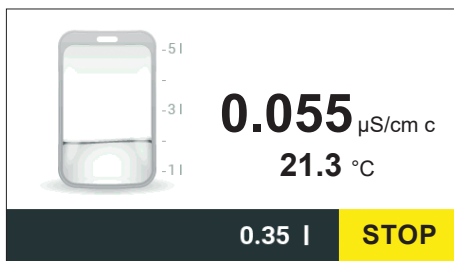
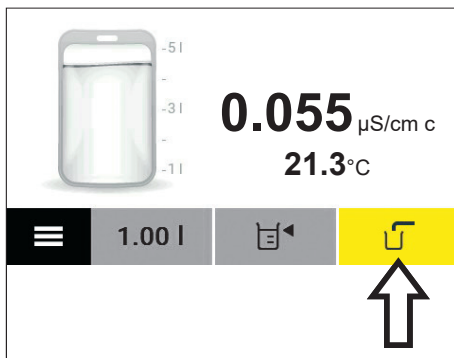
7.2.2 Dispensing Ultrapure Water Manually

With manual dispensing, the ultrapure water is dispensed until the dispensing is stopped manually.

Only Arium® Mini Plus with or without dispensing unit and Arium® Mini: If the fill level of the bag is depleted: Dispensing will end automatically.

Procedure

- ▶ To start the dispensing: Press [Manual dispensing].
 - ▶ Alternatively for devices with a dispensing unit: Press the [Start | Stop] button on the hand-held part.
- ▶ The dispensing starts. The volumes dispensed will be displayed in increments of 0.05 liters (50 ml).
- ▶ The ultrapure water runs into the container with a maximum flow speed of approx. 1.0 l/min.
- ▶ If an Arium® Mini Plus with or without dispensing unit or Arium® Mini is used: The fill level of the bag will be updated on the display during the dispensing.



- ▶ To stop the dispensing: Press the [STOP] button.
 - ▶ Alternatively, for devices with a dispensing unit: Press the [Start | Stop] button on the hand-held part.
- ▶ When the dispensing is complete: Attach the protective cap to the bell assembly of the final filter.

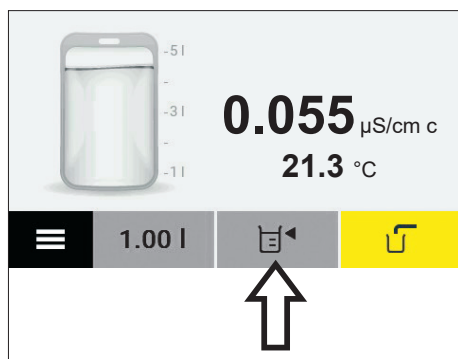
7.2.3 Dispensing Ultrapure Water with Volume-control

During volume-controlled dispensing, a previously specified amount of water is dispensed.

Entering Volumes Using the Numeric Keypad

Procedure

- ▶ Press the [Volume input] symbol.



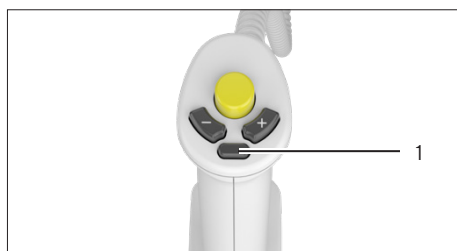
Volume			START
1	2	3	←
4	5	6	
7	8	9	C
0		-	

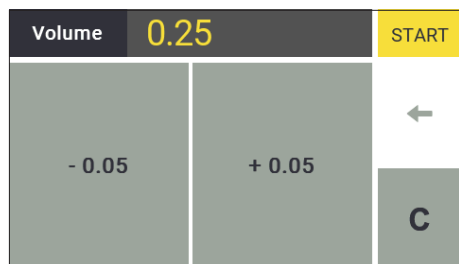
- ▷ The numeric keypad for the volume input appears.
- ▶ Enter the desired dispensing volume in liters. When doing so, observe the requirements for the volume input:
 - The minimum dispense volume is 0.05 liters (50 ml).
 - The maximum dispense volume is 5.00 liters.
 - Only volumes between 0.05 and 5.00 liters can be entered. Invalid numeric fields become inactive (white background).
 - Only Arium® Mini Plus with or without dispensing unit or Arium® Mini: If a higher dispense volume is entered than is available in the bag, the dispensing **cannot** be started. A message to this effect will be sent.
- ▷ The entered volume appears.

Selecting volumes via the hand-held part (for devices with a dispensing unit)

Procedure

- ▶ Press the [Volume] button (1) of the hand-held part.
- ▷ The dialog for the volume selection is displayed.
- ▷ The most recently selected dispensing volume is indicated in the display.





- ▶ To increase the volume: Press the [Plus] button.
- ▶ To decrease the volume: Press the [Minus] button.
- ▷ The volume changes by the amount displayed.
- ▶ If the [Plus] or [Minus] button is pressed briefly: The volume changes by 0.05 liters.
- ▶ If the [Plus] or [Minus] button is pressed and held: The volume changes by 0.5 liters.

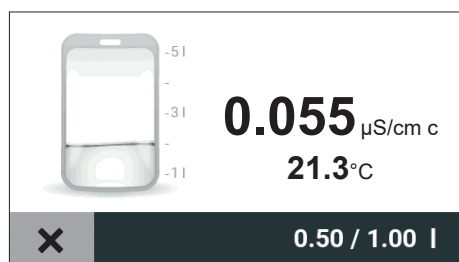
Starting Volume-controlled Dispensing

Procedure

- ▶ Press the [START] button.
 - ▶ Alternatively: Press the [Start | Stop] button on the hand-held part.
- ▷ The dispensing starts:
 - The volumes dispensed will be displayed in increments of 0.05 liters (50 ml).
 - The ultrapure water runs into the container with a maximum flow speed of approx. 1.0 l/min.
 - If an Arium® Mini Plus with dispensing unit or Arium® Mini is used: The fill level of the bag will be updated on the display during the dispensing.
 - The dispensing stops automatically as soon as the selected dispensing volume has been reached.

Canceling Dispensing Prematurely

- ▶ To cancel the dispensing prematurely: Press the [Cancel] button.
 - ▶ Alternatively: Press the [Start | Stop] button on the hand-held part.

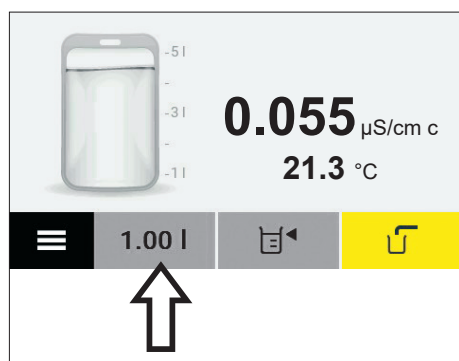


Using Volume-Controlled Dispensing

The last dispensing volumes selected will be saved automatically and displayed on the [Volume-controlled dispense] button on the dispense screen.

Procedure

- ▶ To start another dispensing process with the last dispensing volume selected: Press the [Volume-controlled dispense] button.
- ▷ The dispensing starts.



7.3 Dispensing Ultrapure Water From the Bag (Only With Arium® Mini Plus or Arium® Mini)

If the bag outlet tube is connected to the Bag outlet, ultrapure water can be manually dispensed directly from the bag. Dispensing is carried out exclusively in a depressurized state.

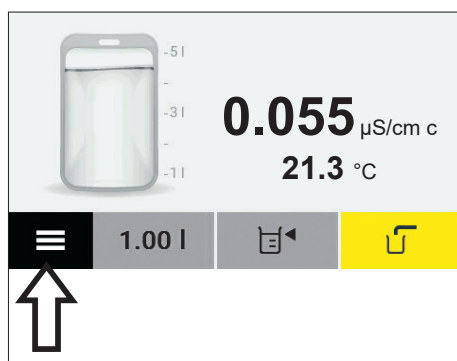
Procedure

- ▶ Place the bag outlet tube into a suitable container.
- ▶ Open the ball valve on the bag outlet tube.
- ▶ The ultrapure water runs out of the bag.
- ▶ When the dispensing is complete: Close the ball valve.

7.4 Opening the Menu

Procedure

- ▶ To open the menu: In dispensing mode, press the [Menu] button. The following tasks can be carried out, e.g. manual bag filling (only Arium® Mini) or opening the “Care” menu.



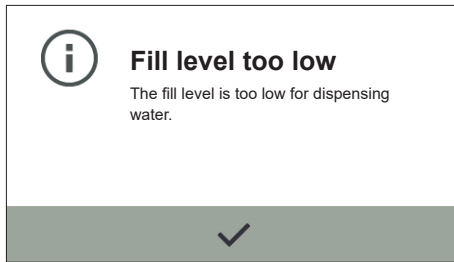
7.5 Manually Fill the Bag (Only for Arium® Mini)

If the bag fill level of the device is depleted or is **no longer** sufficient for the desired dispense volume, the bag can be filled manually. There are various ways in which to begin manual filling of the bag.

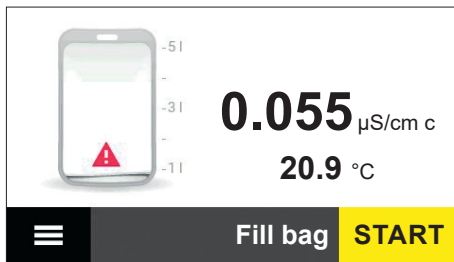
7.5.1 Starting Manual Filling via the Wizard

Procedure

- ▶ If the fill level is too low:
 - ▷ The message “Fill level too low” appears.
 - ▶ Confirm the message with the [Confirmation] button.
 - ▷ The [Fill bag] display appears on the dispense screen.
 - ▶ Press the [START] button.



- ▶ If the bag is already empty prior to dispensing, e.g. due to manual dispensing of ultrapure water:
 - ▷ The [Fill bag] display appears on the dispense screen.
 - ▶ Press the [START] button.



7.5.2 Starting Manual Filling in Menu

If the bag contains water: Bag filling must be launched via the menu. To do so, proceed as follows:

NOTICE

Air ingress may interfere with device functions!

If there is **not** enough water available during the filling process: Air may be drawn in. Air in the system can interfere with the functioning of the device.

- ▶ Prepare a vessel containing a sufficient quantity of pretreated water (at least 5 liters).

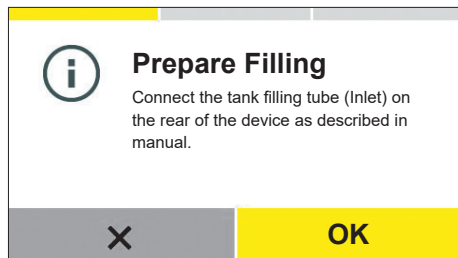
Procedure



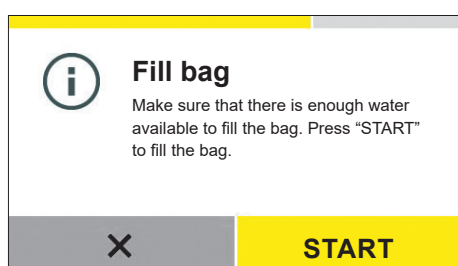
- ▶ Press the [Menu] button.



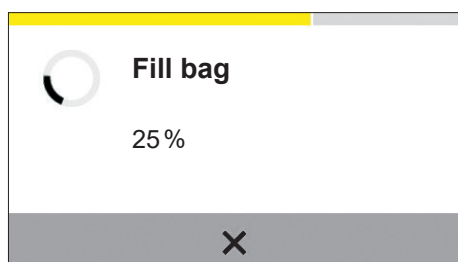
- ▶ Press the [Fill bag] button.



- ▷ The "Prepare filling" (step 1) dialog box appears.
- ▶ Check whether the tank filling tube is connected to the "Inlet" connection.
 - ▶ If required: Connect the tank filling tube.
- ▶ **NOTICE** Contaminated tank filling tube may cause insufficient water quality! Check the tank filling tube for contamination.
 - ▶ If required: Replace the tank filling tube and the bag.
- ▶ Press the [OK] button.



- ▷ The "Fill bag" (step 2) dialog box appears.
- ▶ Place the tank filling tube into the container and secure it to ensure that it is fully submerged throughout the filling process.
- ▶ Press the [START] button.



- ▷ The bag is filled with water. The progress is displayed as a percentage.
- ▶ **NOTICE** Air ingress may interfere with device functions! If there is **not** enough water available, cancel the filling process early and fill the vessel with pretreated water.
- ▶ To cancel the filling process early:
 - ▶ Press the [Cancel] button.
- ▷ The "Fill bag" (step 2) dialog box appears.
- ▶ Carry out the filling procedure again as described above.
- ▷ When the filling procedure is complete, the "Filling finished" dialog box appears.
- ▶ Check whether the free end of the tank filling tube is stored securely.
 - ▶ If required: Disconnect the tank filling tube from the "Inlet" connection on the rear of the device and store in a safe place.
- ▶ Confirm the filling process by pressing the [OK] button.
- ▷ The display changes to the dispensing screen.

7.6 Activating or Deactivating Standby Mode

If **no** water has been dispensed for an extended period, the device automatically switches to standby mode. This ensures economical and eco-friendly operation. In standby mode, the pretreatment stage is **no longer** active and the water in the ultrapure water stage circulates in intervals. Standby mode can also be activated manually.

If the device is manually switched to standby mode while the bag is being filled automatically, the user will be asked if they wish to proceed. The bag filling can be canceled directly, putting the device into standby mode. If the filling is continued, the device automatically switches to standby mode afterwards.

Procedure



- ▶ To activate standby mode: Press the [Standby] button.
- ▷ The display goes dark. The backlight on the [Standby] button stays visible.
- ▶ To deactivate standby mode: Press the [Standby] button.
- ▷ The start screen appears. Once the system start is complete, the display changes to the dispensing screen.

7.6.1 Automatic ECO Mode

As well as standby mode, the device also has an automatic ECO mode. One minute after the final procedure, recirculation of the ultrapure water stops and the display darkens. After a further 15 minutes without anyone touching the display, the device automatically switches to standby mode.

7.7 Changing System Settings

The system settings for the device, e.g., date, time, displayed values, can be changed in the “Settings” menu.

Procedure



- ▶ Press the [Menu] button.



- ▶ Press the [Settings] button.
- ▷ The “Settings” menu appears.
- ▶ Configure the desired settings (for possible settings, see the parameter list (see Chapter “4.10 Parameters of the “Settings” Menu,” page 28).

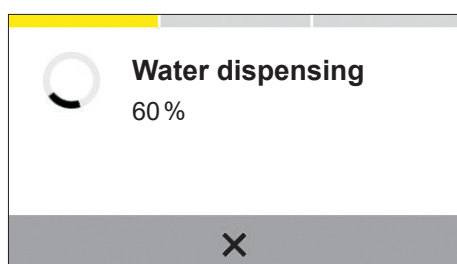
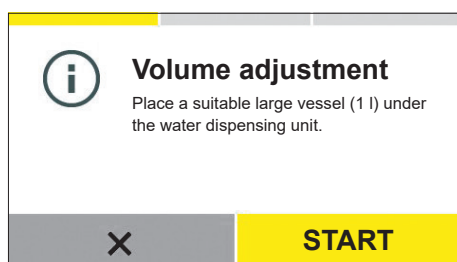
7.8 Performing a Volume Adjustment

The flow sensor of the device can be readjusted to be able to dispense the most precise possible volume of ultrapure water.

A sample volume of around half a liter is dispensed for this purpose. The actual volume of the dispensed sample volume is measured and transmitted to the device. This sample volume is used as a reference for the volume dispensing.

Procedure

- ▶ Press the [Volume adjustment] menu option.
- ▶ The “Volume adjustment” dialog box appears.
- ▶ Place a graduated measuring beaker or cylinder with a capacity of at least 1 liter under the water outlet.
- ▶ Press the [START] button.



- ▶ The “Water dispensing” dialog box appears. The output progress is displayed as a percentage.
- ▶ Once 100 % has been reached, the dispensing process stops automatically.
- ▶ The “Water volume” dialog box appears. The liter value of the sample volume (output water volume) must be entered.

Using the Weight of the Sample Volume to Determine the Volume

If a graduated measuring beaker or cylinder is **not** available, the weight of the sample volume can be used as an alternative to determine the actual volume.

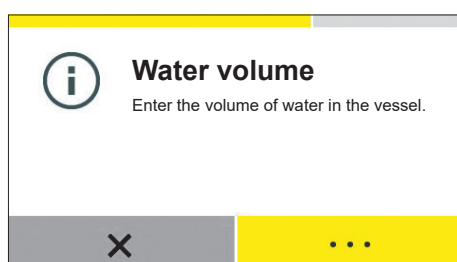
Procedure

- ▶ Determine the weight of the sample volume.
- ▶ Convert the weight of the sample volume into its value in liters (1 g ultrapure water is equivalent to 1 ml ultrapure water).

Enter the Sample Volume in Liters

Procedure

- ▶ Press the [Input] button.
- ▶ Enter the liter value of the sample volume.
- ▶ The “Adjustment finished” dialog box appears.
- ▶ Confirm the procedure with the [Confirmation] button.



8 Cleaning and Maintenance

8.1 Cleaning

8.1.1 Cleaning the Display

Procedure

- ▶ To **avoid** uncontrolled changes to the settings of the device:
Activate standby mode (see Chapter 7.6, page 58).
- ▶ Wipe the display gently with a soft, dry cloth.
- ▶ Deactivate standby mode (see Chapter 7.6, page 58).

8.1.2 Cleaning the Device Housing

CAUTION

Danger of injury caused by electric current!

When cleaning the device and its components while they are connected to the power supply, there is a risk of electric shock.

- ▶ Always disconnect the device from the AC power before cleaning.

NOTICE

The electronic equipment could be damaged by improper cleaning!

Liquids or dust can damage the device or the power supply.

- ▶ **Never** open the power supply or the right side of the device housing.
- ▶ Make sure that **no** liquids or dust get into the device or the AC adapter.

NOTICE

Damage to the device surfaces!

Aggressive cleaning agents may damage device surfaces.

- ▶ Never use cleaning agents that contain solvents, acetone or abrasive ingredients.

Procedure

- ▶ Disconnect the device from the power supply.
- ▶ Wipe the housing of the device with a damp cloth.
- ▶ If installed components must be cleaned: Remove the front cover and the left side cover and wipe down the installed components with a damp cloth.
- ▶ Dry the device with a soft cloth.

8.2 Maintenance Schedule

Depending on the volume of water dispensed, it may be necessary to change the consumables more often than specified in the maintenance schedule. If, for example, sterile water is required, the final filter must be replaced regularly.

It is advisable to replace different consumables in the same maintenance step. This saves time and water. We recommend that you make a sensible plan for replacing consumables.

Interval	Component	Task	Chapter, page
12 months	UV lamp (optional)	Changing the UV Bulb	8.5.5, page 66
Max. 6 months (depending on the volume of water dispensed)	Ultrapure water cartridge (Scientific Pack)	Replace ultrapure water cartridge	8.5.4, page 65
6 months (only Arium® Mini Plus with or without dispensing unit or Arium® Mini)	Bag	Replacing the Bag	8.5.2, page 63
6 months (only Arium® Mini Plus with or without dispensing unit)	Pretreatment cartridge	Replace the pretreatment cartridge	8.5.3, page 64
1 – 24 weeks (depending on the application)	Sterile filter (finalfilter)	Change final filter	8.5.6, page 69
1 – 13 weeks (depending on the application)	Ultrafilter (final filter)	Change final filter	8.5.6, page 69

8.3 Opening the “Care” Menu

The “Care” menu contains all the steps for the care and maintenance work.

Procedure



- ▶ Press the [Care] symbol in the menu.
- ▷ The “Care” menu appears, and tasks such as the following can be carried out: Display reminders, replace consumables.

8.4 Displaying Reminders

Reminders to replace certain consumables are automatically displayed as warning messages (see Chapter 9.2, page 76). Pending consumables replacements can be viewed at a glance.

All reminders are automatically updated after the consumable has been replaced.

Procedure

- ▶ Tap the menu item [Reminder] in the “Care” menu.
- ▷ The date of the next required replacement is displayed, e.g. for the bag, the cleaning cartridge or the UV lamp.

Reminder		✓
Bag:	01.07.2016	
Pretreatment cartridge:	01.07.2016	
Ultrapure water cartridge:	01.07.2016	
UV lamp:	01.02.2017	
Sterile final filter:	01.04.2016	

8.5 Replacing Consumables

8.5.1 Selecting Consumables for Replacement

Consumable replacement must be set using the display. To do this, you must define which consumables are to be replaced in the menu “Replace consumable”. One or more consumables can be selected.

If several consumables are selected: The wizard guides you through the replacement of the individual consumables. The individual steps are shown on the display.

Procedure

- ▶ Tap the menu item [Change material] in the “Care” menu.
- ▷ The “Change material” menu appears. The consumables that can be changed are indicated, e.g. bag, pretreatment cartridge.
- ▶ Select the desired consumables. To do this, tap on the menu items for the consumables, e.g. [Bag], [Pretreatment cartridge] and [Ultrapure water cartridge].
- ▷ The selected menu items are highlighted and marked with a checkmark.
- ▶ Confirm the selection of the consumables with the [Confirmation] button.
- ▷ If several consumables are selected: The wizard guides you through all consumable replacement and work steps.

Change material		✓
✓	Bag	
✓	Pretreatment cartridge	
✓	Ultrapure water cartridge	
	UV lamp	
	Sterile filter	

8.5.2 Replacing Bag (only Arium® Mini Plus or Arium® Mini)

Draining the Bag

Requirements

- In the “Change material” menu, the menu item [Bag] is activated.
- The “Empty bag” dialog box appears.

Procedure

- ▶ Check whether the tank outlet tubing is connected to the “Bag Outlet” connection.
 - ▶ If required: Connect the tank outlet tubing.
- ▶ Guide the tank outlet tubing into a suitable container.
- ▶ In order to empty the bag completely: Place the container underneath the device.
- ▶ Open the ball cock of the tank outlet tubing.
- ▷ Pure water flows out of the bag.
- ▶ When the bag is empty: Confirm the emptying of the bag with the [OK] button.
- ▷ The dialog box “Change bag” appears.

Removing Empty Bag

Procedure

- ▶ Remove the left side cover (see Chapter “5.3.1 Removing the Left Side Cover,” page 32).
- ▶ To make it easier to reach the lower connections on the device: Pull the tray up and remove it from the device.
- ▶ Disconnect the gray quick connector on the three device connections successively.
- ▶ Slide the empty bag out of the guide rail and remove it from the device.



Inserting New Bag

- ▶ Insert a new bag (see Chapter “6.5 Inserting Bag (only Arium® Mini Plus or Arium® Mini),” page 41).
- ▷ The dialog box “Close tank outlet” appears.
- ▶ Close the ball cock of the tank outlet tubing.
- ▶ Confirm that the tank outlet has been closed with the [OK] button.

- ▷ If an Arium® Mini Plus with or without dispensing unit is used: The device automatically begins filling the bag.
- ▶ If an Arium® Mini is used: The filling process must be started.
 - ▷ The dialog box “Fill bag?” appears on the device display.
 - ▶ To fill the bag: Tap the [YES] button. Manually fill the bag (see Chapter 7.5.2, page 56).
 - ▶ To fill the bag at a later point in time: Tap the [NO] button.

8.5.3 Replacing Pretreatment Cartridge (only Arium® Mini Plus)

To replace pretreatment cartridges, the final filter must be removed and the device depressurized. The pretreatment cartridge is marked with an “R”.

Requirements

- In the “Change material” menu, the menu item [Pretreatment cartridge] is activated.
- The dialog box “Remove filter” appears.

Procedure

- ▶ Remove the final filter (removal, see Chapter “8.5.6 Changing Final Filter,” page 69).
- ▷ The dialog box “Prepare depress. / Start depressurization” appears.
- ▶ Start depressurization (see Chapter 8.7, page 71).
- ▷ When depressurization is complete, the dialog box “Remove cover” appears.
- ▶ Remove the front cover.
- ▶ Confirm removal of the cover with the [OK] button.
- ▷ The dialog box “Change cartridge” appears. The pretreatment cartridge (R) can be changed.
- ▶ Press the two protruding clamps of the pretreatment cartridge together and pull the pretreatment cartridge forward and out.



- ▶ Confirm removal of the pretreatment cartridge with the [OK] button.
- ▶ Insert a new pretreatment cartridge in the device (see Chapter 6.3, page 39).
- ▶ Rinse the pretreatment cartridge (see Chapter 6.7, page 45).
- ▷ Once the pretreatment cartridge has been rinsed: The dialog box “Connect filter” appears.
- ▶ Connect the final filter (see Chapter 6.9, page 49).
- ▷ Replacing the pretreatment cartridge is complete.

8.5.4 Replacing Ultrapure Water Cartridge

To replace the ultrapure water cartridge, the final filter must be removed and the device depressurized. The ultrapure water cartridge is marked with an “L”.

Requirements

- In the “Change consumable” menu, the menu item [Ultrapure water cartridge] is activated.
- The dialog box “Remove filter” appears.

Procedure

- ▶ Remove the final filter (removal, see Chapter “8.5.6 Changing Final Filter,” page 69).
- ▷ The dialog box “Prepare depress. / Start depressurization” appears.
- ▶ Start depressurization (see Chapter 8.7, page 71).
- ▷ When depressurization is complete, the dialog box “Remove cover” appears.
- ▶ Remove the front cover.
- ▶ Confirm removal of the cover with the [OK] button.
- ▷ The dialog box “Change cartridge” appears. The ultrapure water cartridge (L) can be changed.
- ▶ Press the two protruding terminals of the ultrapure water cartridge together and pull the ultrapure water cartridge forward and out.



- ▶ Confirm removal of the ultrapure water cartridge with the [OK] button.
- ▶ Insert a new ultrapure water cartridge in the device (see Chapter 6.4, page 40).
- ▶ Rinse the ultrapure water cartridge (see Chapter 6.8, page 46).
- ▷ Once the ultrapure water cartridge has been rinsed: The dialog box "Connect filter" appears.
- ▶ Connect the final filter (see Chapter 6.9, page 49).
- ▷ Replacing the ultrapure water cartridge is complete.

8.5.5 Replacing the UV Lamp

The device can be equipped with a UV lamp. When replacing the UV lamp: A suitable UV lamp must be used (suitability, see Chapter "15 Consumables," page 89).

The UV lamp is installed under the right-side cartridge holder (pretreatment cartridge). The pretreatment cartridge is marked with an "R".

To change the UV lamp, the final filter must be removed and the device depressurized.

If an Arium® Mini Plus with or without dispensing unit is used: The pretreatment cartridge must be removed to access the UV lamp behind it.

Requirements

- In the "Change material" menu, the menu item [UV lamp] is activated.
- The dialog box "Remove filter" appears.

Procedure

- ▶ Remove the final filter (removal, see Chapter "8.5.6 Changing Final Filter," page 69).
- ▷ The dialog box "Prepare depress. / Start depressurization" appears.
- ▶ Start depressurization (see Chapter 8.7, page 71).
- ▷ When depressurization is complete, the dialog box "Remove cover" appears.
- ▶ Remove the front cover.
- ▶ Confirm removal of the cover with the [OK] button.
- ▶ If an Arium® Mini Plus with or without dispensing unit is used:
 - ▷ The "Remove the pretreatment cartridge (R) as described in the manual" dialog box appears.
 - ▶ Remove the pretreatment cartridge (see Chapter 8.5.3, page 64).
- ▷ The dialog box "Change UV lamp" appears.

Removing the UV Lamp

⚠ CAUTION

Risk of injury from electrical current and UV radiation!

The UV lamp emits UV radiation and may be live.

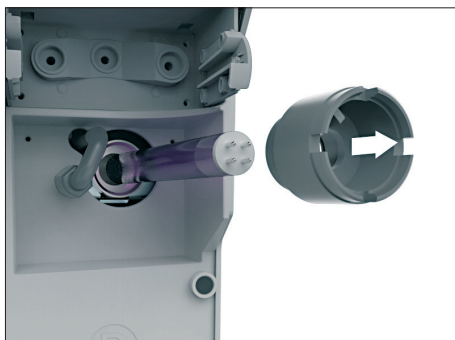
- ▶ Disconnect the device from the power supply before the old UV lamp is removed.

Procedure

- ▶ Disconnect the device from the power supply.
- ▶ The display goes out. After the power has been restored, the wizard will continue automatically.
- ▶ Squeeze the metal retaining clip on the old UV lamp and pull it forward to remove it.
- ▶ Remove the metal retaining clip from the cable and store it in a safe place, e.g. on the magnet securing the cover.



- ▶ Disconnect the black connector from the old UV lamp.
- ▶ Unscrew the black plastic cover of the UV lamp housing. If required: Use a suitable tool to help you do this.



- ▶ Remove the black plastic cover of the UV lamp housing.
- ▶ Carefully pull the old UV lamp out of the lamp unit and dispose of it properly (see Chapter "13 Disposal," page 81).

Inserting the UV Lamp

NOTICE

Touching UV lamps with your bare fingers will cause them to become defective!

Touching the UV lamp with your bare fingers will leave fingerprints. The fingerprints can become so hot during operation that the UV lamp is destroyed.

- ▶ **Never** touch the glass of the UV lamp with your bare fingers.
 - ▶ Only hold the UV lamp where it connects to the device or when wearing gloves.
-

Procedure

- ▶ Carefully unpack the new UV lamp without touching the glass with your fingers.
- ▶ Insert the new UV lamp all the way into the lamp housing without exerting pressure on it.
- ▶ Unscrew the black plastic cover of the UV lamp housing by hand.
- ▶ Attach the black connector to the UV lamp. The connector only fits in two orientations. Both orientations allow the UV lamp to function.
- ▶ Slide any protruding cable back into the housing.
- ▶ Slide the metal retaining clip over the cable and onto the black plastic cover of the UV lamp housing.
- ▶ Insert the closed side of the metal retaining clip into the recess of the black plastic cover provided.
- ▶ Squeeze the open side of the metal retaining clip and insert it into the recesses provided in the black plastic cover.
- ▷ Release the metal retaining clip secures it to the black plastic cover.
- ▶ Connect the device to the power supply.
- ▷ The wizard continues.
- ▶ Follow the instructions on the display, e.g. "Insert the new pretreatment cartridge (R) as described in the manual".

8.5.6 Changing Final Filter

Remove final filter

Requirements

- If the [Sterile filter] or [Ultrafilter] menu item is activated in the “Replace consumables” menu: The “Remove filter” dialog box appears.
- If the menu items [pretreatment cartridge], [ultrapure water cartridge] or [UV lamp] are activated in the “Replace consumable” menu: The dialog box “Remove filter” appears.

Procedure



- ▶ Use the tubing removal tool to push and hold the water outlet quick connector up.
- ▷ The quick connector is unlocked.



- ▶ For devices with a dispensing unit: Use the tubing release tool to push up and hold the quick connector of the water outlet on the hand-held part.
- ▷ The quick connector is unlocked.
- ▶ Pull the final filter out of the quick connector.
- ▶ Confirm removal of the final filter with the [OK] button.

Connect new final filter

Requirements

- The final filter has been removed.
- The dialog box “Connect filter” appears.

Procedure

- ▶ Connect and rinse a new final filter (see Chapter 6.9, page 49 and Chapter 6.10, page 50).

8.6 Enabling, Disabling or Configuring Reminders for Replacing Final Filters

If sterile or endotoxin-free water is constantly required, the final filter must be replaced regularly. The device can provide a reminder about a pending final filter replacement.

Procedure

- ▶ Tap the menu item [Final filter reminder] in the “Care” menu.
- ▶ The dialog box “Final filter reminder” appears.
- ▶ Tap the [Active] button.
- ▶ Select the installed final filter, e.g. [sterile filter] or [ultrafilter].
- ▶ The currently configured replacement interval appears in the column “Reminder [weeks]”.
- ▶ To change the replacement interval: Press the [Enter] button.
- ▶ The numeric keypad appears.
- ▶ Type in the desired replacement interval (in weeks) (for required maintenance intervals, see Chapter “8.2 Maintenance Schedule,” page 61).
- ▶ Confirm the input with the [Confirm] button.
- ▶ To activate the reminder: In the “Final filter reminder” dialog box, press the [Confirmation] button.
- ▶ If **no** sterile or endotoxin-free water is required:
 - ▶ Remove the final filter (see Chapter 8.5.3, page 64).
 - ▶ To deactivate reminders about replacing the final filter: In the dialog box “Final filter reminder”, press the [Inactive] button.

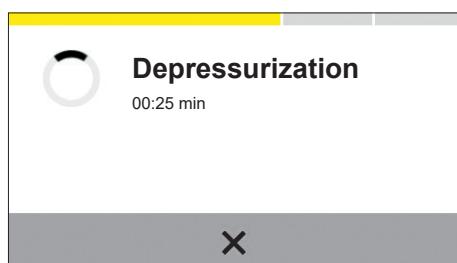
Final filter reminder		
Final filter reminder	Filter Type	Reminder [weeks]
Active	Sterile filter	4
Inactive	Ultrafilter	...

8.7 Carrying Out Depressurization

The device is under pressure during operation. If the device is taken out of service for an extended period of time or permanently, the pressure in the device must be let out manually.

Procedure

- ▶ Tap the menu item [Depressurize] in the “Care” menu.
- ▷ The dialog box “Remove filter” appears.
- ▶ Remove final filter (see Chapter “8.5.6 Changing Final Filter,” page 69).
- ▶ Confirm removal of the final filter with the [OK] button.
- ▷ The dialog box “Prepare depress.” appears.
- ▶ Place a vessel (at least 1 liter) under the water outlet. Alternatively, connect the dispense tube and run the free end to the drain.
- ▶ Confirm this preparatory step with the [OK] button.
- ▷ That dialog box “Start depressurization” appears.
- ▶ Tap the [START] button.
- ▷ The dialog box “Depressurization” appears.
- ▷ The device depressurizes. The process takes about half a minute.
- ▶ To cancel the depressurization process before it is finished, e.g., if there is **not** a sufficiently large vessel available: Press the [Cancel] button.
- ▷ That dialog box “Start depressurization” appears again.
- ▶ Carry out depressurization as described above again.
- ▷ When depressurization is complete, the dialog box “Turn off device” appears.
- ▶ Disconnect the device from the power supply.
- ▶ Pressurization is carried out when the device is switched on again.
- ▶ Perform the rinsing process (see Chapter “8.9 Performing the Final Rinsing Procedure,” page 72)



8.8 Carrying Out Venting

During venting, the device fills and rinses the ultrapure water cartridge. Air is removed from the ultrapure water circulation in the process.

Given the following conditions: The ultrapure water circuit must be purged:

- The displayed water quality is fluctuating continuously during operation.
- The device was taken out of service for an extended period.

Procedure

- ▶ Tap the menu item [Venting] in the “Care” menu.
- ▷ The dialog box “Remove filter” appears.
- ▶ Remove the final filter (see Chapter 8.5.3, page 64).
- ▶ Confirm removal of the final filter with the [OK] button.
- ▷ The dialog box “Prepare rinsing” appears.
- ▶ Place a vessel (at least 1 liter) under the water outlet. Alternatively, connect the dispense tube and run the free end to the drain.
- ▶ Tap the [OK] button.

- ▶ Arium® Mini Plus with or without dispensing unit product water quality:
 - ▷ If the bag **no** longer contains a sufficient amount of water or empties during the rinsing process, the bag is filled automatically.
 - ▶ When filling is complete, the dialog box “Preparing to rinse” appears: Start the rinsing process again.
- ▶ Only Arium® Mini:
 - ▷ If the bag **no** longer contains a sufficient amount of water or empties during the rinsing process, the dialog box “Fill bag” appears.
 - ▶ The dialog box “Fill bag” appears: Manually fill the bag (see Chapter 7.5.2, page 56).

- ▶ Perform the rinsing process (see Chapter “8.9 Performing the Final Rinsing Procedure,” page 72).

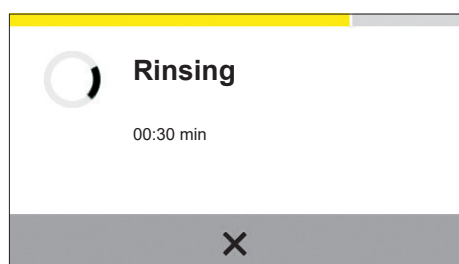
8.9 Performing the Final Rinsing Procedure

After depressurizing or venting, the device needs to be rinsed.

Procedure

- ▷ The “Prepare rinsing” dialog box appears.
- ▶ Press the [OK] button.
- ▷ The “Start rinsing” dialog box appears.
- ▶ Press the [START] button.
- ▷ During the rinsing process, the remaining rinse time is displayed in minutes.
- ▶ To interrupt the rinsing process:
 - ▶ Press the [Cancel] button.
 - ▷ The “Start rinsing” dialog box is displayed again.
- ▶ To resume the rinsing process after an interruption: Press the [START] button.

- ▷ When rinsing is complete, the “Connect filter” dialog box appears.
- ▶ Connect the final filter (see Chapter 6.9, page 49).



9 Malfunctions

9.1 Error Messages

If an error message is active, dispensing is canceled and locked automatically.

Error message	Fault	Cause	Correction	Chapter, page
Error 0105	Measured values are not being displayed.	There is an internal communications error.	Disconnect the device from the AC power and wait a minute. Reconnect the device to the AC power. If the error persists: Contact Sartorius Service.	7.1, page 51
Error 0140	The UV lamp is not recognized.	The UV lamp is not connected properly or defective.	Check whether the black plug of the UV lamp is connected correctly. Remove the UV lamp and check for damage. If required: Change the UV lamp. If the error persists: Contact Sartorius Service.	8.5.5, page 66
Check UV lamp, replacement may be needed.				8.5.5, page 66
Error 0150	The device has a leak.	Water has leaked inside the device.	Remove the left side cover and check that the three bag hoses are properly connected to the device connectors. If required: Remove the bag and reconnect it. Drain the tray at the bottom of the device and check that the connections are tight during operation. If the error persists: Contact Sartorius Service.	5.3.1, page 32
Check device for leakage. Please contact your service technician.				
Error 0160	Operation of the device is not possible.	The front cover is not properly attached.	Correctly attach the front cover to the device. If the error persists: Contact Sartorius Service.	
Front cover is open. Please put the front cover back on.				

Error message	Fault	Cause	Correction	Chapter, page
Error 0163 Check proper fit of the pretreatment cartridge (R).	The pretreatment cartridge is not recognized.	The pretreatment cartridge has not been properly inserted.	Press the pretreatment cartridge in firmly until you hear a distinct clicking sound. If the error persists: – Remove the pretreatment cartridge. – Insert the pretreatment cartridge.	8.5.3, page 64
Error 0166 Check proper fit of the ultrapure water cartridge (L).	The ultrapure water cartridge is not detected.	The ultrapure water cartridge has not been properly inserted.	Press the ultrapure water cartridge in firmly until you hear a distinct clicking sound. If the error persists: – Remove the ultrapure water cartridge. – Insert the ultrapure water cartridge..	8.5.4, page 65
Error 0180 Please contact your service technician.	The fill level of the bag is not being properly detected.	The bag is damaged or improperly connected. The sensor is defective.	Remove the left side cover and check the bag for damage. If required: Change bag. Check that the three bag hoses are properly connected to the device connectors. If required: Remove the bag and reconnect it. Using the tank outlet tubing connected to "Bag Outlet", let about ¼ liters of pure water out of the bag and check the level change.	5.3.1, page 32 8.5.2, page 63 7.3, page 55
Error 0206	No connection to the hand-held part.	– The device cannot establish a connection to the hand-held part. – The hand-held part connection cable is not connected.	Connect the connection cable for the dispensing unit to the device.	5.4.1, page 34

Error message	Fault	Cause	Correction	Chapter, page
Error The conductivity of the ultrapure water is outside of the measuring range.	The conductivity of the ultrapure water is outside of the measuring range.	There is air in the ultrapure water circulation.	Initiate venting.	8.8, page 71
		The ultrapure water cartridge has been used up.	Replace the ultrapure water cartridge.	8.5.4, page 65
			To dispense water for testing purposes: Disable the dispensing lock. If the error persists: Contact Sartorius Service.	
Error The temperature of the ultrapure water is outside of measuring range.	The temperature of the ultrapure water is outside of measuring range.	The temperature of the ultrapure water is outside of measuring range.	Check whether the ambient temperature complies with the device specifications in the technical data.	14.2, page 83
			Dispense about 1 liter: – To dispense water for testing purposes: Disable the dispensing lock. – Dispense and discard about 1 liter of water.	
			If the error persists: Contact Sartorius Service.	
Error Ultrapure water quality > XX µS/cm	The limit has been exceeded and the dispensing lock is preventing further dispensing.	There is air in the ultrapure water circulation.	Perform venting.	8.8, page 71
			Replace the ultrapure water cartridge.	8.5.4, page 65
			To dispense water for testing purposes: Disable the dispensing lock. If the error persists: Contact Sartorius Service.	
		The limit has been configured incorrectly.	Check the limit. If required: Configure the limit.	

9.2 Warning Messages

If a warning message is active, water can still be dispensed. The water quality is impaired under certain circumstances.

Warning message	Fault	Cause	Correction	Chapter, page
The conductivity measurement of the RO water is out of range.	The conductivity of the pure water cannot be determined.	The quality of the feed water is insufficient .	Check whether the quality of the feed water complies with the device specifications in the technical data.	14.6, page 86
		The pretreatment cartridge has been used up.	Replace the pretreatment cartridge. If the error persists: Contact Sartorius Service.	8.5.3, page 64
The conductivity of the ultrapure water is outside of the measuring range.	The conductivity of the ultrapure water cannot be determined.	There is air in the ultrapure water circulation.	Perform venting.	8.8, page 71
		The ultrapure water cartridge has been used up.	Replace the ultrapure water cartridge. If the error persists: Contact Sartorius Service.	8.5.4, page 65
Temperature measurement of the RO water is out of range.	The temperature of the pure water cannot be determined.	The temperature of the pure water is outside of measuring range.	Check whether the temperature of the feed water complies with the device specifications in the technical data.	14.6, page 86
			Check whether the ambient temperature complies with the device specifications in the technical data.	14.2, page 83
			If the error persists: Contact Sartorius Service.	
The temperature of the ultrapure water is outside of measuring range.	The temperature of the ultrapure water cannot be determined.	The temperature of the ultrapure water is outside of measuring range.	Check whether the ambient temperature complies with the device specifications in the technical data.	14.2, page 83
			Dispense and discard about 1 liter of water.	
			If the error persists: Contact Sartorius Service.	
RO water quality > XX µS/cm	The limit has been exceeded.	The quality of the pure water is insufficient .	Check whether the quality of the feed water complies with the device specifications in the technical data.	14.6, page 86
		The pretreatment cartridge has been used up.	Replace the pretreatment cartridge.	8.5.3, page 64

Warning message	Fault	Cause	Correction	Chapter, page
Ultrapure water quality > XX µS/cm	The limit has been exceeded.	There is air in the ultrapure water circulation.	Perform venting.	8.8, page 71
			Replace the ultrapure water cartridge.	8.5.4, page 65
		The limit has been configured incorrectly.	Check the limit. If required: Configure the limit.	
RO water temperature > XX °C	The limit has been exceeded.	The temperature of the feed water is too high or too low.	Check whether the temperature of the feed water complies with the device specifications in the technical data.	14.6, page 86
Ultrapure water temperature > XX °C	The limit has been exceeded.	The ambient temperature of the device is too high or too low.	Check whether the ambient temperature complies with the device specifications in the technical data.	14.2, page 83
			Dispense and discard about 1 liter of water.	
Replacement of the pretreatment cartridge (R) is required.	The pretreatment cartridge needs to be replaced.	The replacement interval of the pretreatment cartridge has expired.	Replace the pretreatment cartridge.	8.5.3, page 64
Replacement of the ultrapure water cartridge (L) is required.	The ultrapure water cartridge needs to be replaced.	The replacement interval of the ultrapure water cartridge has expired.	Replace the ultrapure water cartridge.	8.5.4, page 65
Replacement of the bag is required.	The bag needs to be replaced.	The replacement interval of the bag has expired.	Replace the bag.	8.5.2, page 63
Replacement of the UV lamp is required.	The UV lamp needs to be replaced.	The replacement interval of the UV lamp has expired.	Change UV lamp.	8.5.5, page 66
Replacement of the sterile final filter is required.	The sterile final filter needs to be replaced.	The replacement interval of the sterile final filter has expired.	Replace the final filter.	8.5.3, page 64
Sartorius maintenance is required.	Maintenance service must be performed.	The maintenance service interval has expired.	Contact Sartorius Service.	

9.3 Additional Faults

Fault	Cause	Correction	Chapter, page
The device unexpectedly stops dispensing.	The bag is empty.	Check the level of the bag on the display.	
		Remove the left cover and check the fill level of the bag. If required: Manually fill the bag (only Arium® Mini).	7.4, page 55
		Check that the feed water tubing has been correctly connected.	6.6.1, page 42
	There is no feed water is connected.	Connect the feed water tubing or tank filling tubing.	6.6, page 42
	The final filter is clogged or contains air.	If the final filter is attached and no water can be dispensed:	
		– Rinse the final filter.	6.10, page 50
		– Vent the final filter.	8.8, page 71
		– If the error persists: Replace the final filter.	8.5.6, page 69
		If the final filter has been removed and no water can be dispensed: Contact Sartorius Service.	

10 Storage and Shipping

10.1 Storage

Procedure

- ▶ If the device is in operation:
 - ▶ Decommission the device.
 - ▶ Clean the device.
- ▶ Store the device according to the ambient conditions (see Chapter "14.2 Ambient Conditions," page 83).

10.2 Returning Device and Parts

You can send defective devices or parts back to Sartorius. Returned devices must be clean, decontaminated and properly packed.

Transport damage as well as measures for subsequent cleaning and disinfection of the device or parts by Sartorius shall be charged to sender.

WARNING

Risk of injury due to contaminated equipment!

Devices contaminated with hazardous materials (NBC contamination) will **not** be accepted for repair or disposal.

- ▶ Observe the information on decontamination (see Chapter 13.1, page 81).
-

Procedure

- ▶ Decommission the device.
- ▶ Contact Sartorius Service for instructions on how to return equipment or parts (please refer to our website at www.sartorius.com for return instructions).
- ▶ Pack the device and its parts properly for return.

11 Decommissioning

Requirement

Operation has been ended correctly.

Procedure

- ▶ Start depressurization (see Chapter 8.7, page 71)
- ▶ Disconnect the device from power.
- ▶ Disconnect the device from the supply lines. Remove the consumables being used.
- ▶ Disconnect any attached components from the device.
- ▶ Clean the device (see Chapter 8.1, page 60).

12 Transport

12.1 Transporting the Device

Requirements

The device has been taken out of operation.

Procedure

- ▶ **⚠ CAUTION** Injury to the feet due to falling device components during improper transport!
 - ▶ Do not carry the device using the components of the dispensing unit.
 - ▶ **Do not** lift or transport the device by the side covers.
 - ▶ Hold the device underneath the display on the front side and by the recess for the power supply unit on the back side and lift it carefully.



13 Disposal

13.1 Information on Decontamination

The device does **not** contain any hazardous materials that would necessitate special disposal measures.

If the device has come into contact with hazardous substances: Measures must be carried out to properly decontaminate and declare such devices. The operator is responsible for adhering to local legislation on the proper declaration of transport and disposal and the proper disposal of the device.

WARNING

Risk of injury due to contaminated equipment!

Devices contaminated with hazardous materials (NBC contamination) will **not** be accepted by Sartorius for repair or disposal.

13.2 Disposing of Device and Parts

13.2.1 Information on Disposal

The device and its accessories do **not** belong in regular household waste, since they are made of high-grade materials that can be recycled and reused. All parts must be disposed of properly by disposal facilities.

Batteries are installed in the device. Batteries do **not** belong in regular household waste, since they are made of high-grade materials that can be recycled and reused. Batteries must be disposed of properly by disposal facilities.

The packaging is made of environmentally friendly materials that can be used as secondary raw materials.

The consumables are designed and intended for single-use.

Hazardous Substances

The UV lamp contains mercury. The UV lamp must be delivered to an approved disposal center for hazardous substances.

13.2.2 Disposal

Requirements

If the device has come into contact with hazardous substances: The device and consumables have been decontaminated.

Procedure

- ▶ Remove the UV lamp from the device (see Chapter “8.5.5 Replacing the UV Lamp,” page 66).
- ▶ Deliver the UV lamp to an approved disposal center for hazardous substances.
- ▶ Dispose of the device. Follow the disposal instructions on our website (www.sartorius.com). Inform the disposal facility that batteries are installed in the device.
- ▶ Dispose of the packaging in accordance with local government regulations.
- ▶ Dispose of the consumables in accordance with local government regulations.

14 Technical Specifications

14.1 Power Supply

	Unit	Value
Sartorius power supply, Model 1000018304		
Primary		
Voltage	V~	100 - 240 (±10%)
Frequency	Hz	50 to 60
Current, max.	A	2.0
Secondary		
Voltage	V=	+24 (<5%)
Current, max.	A	6.25
Short circuit protection		Electronic
Protection class		I
Height above sea level, maximum	m	3000
Pollution level of the power supply unit according to IEC 61010-1		2
Power supply connection cable		
Connection plug according to DIN EN/IEC 60320-1/C14		Country-specific, 3-pin, two-sided plug
Connector according to DIN EN/IEC 60320-1/C14		3-pin
Other data		See power supply label
Device		
Power supply		Only via Sartorius power supply 1000018304
Input supply voltage	V _{DC}	+24 (±10%)
Current consumption, max.	A	3.0

14.2 Ambient Conditions

	Unit	Value
Environment		For indoor use only
Storage and shipping temperature	°C	+5 to +45
Operating temperature	°C	+2 to +35
Operating height	m above sea level	Up to 3000
Relative humidity	%	40 to 80

14.3 Safety of Electrical Equipment

Safety requirements according to
DIN EN/IEC 61010-1

Safety requirements for electrical equipment for
measurement, control and laboratory use – Part 1:
General requirements

14.4 Electromagnetic Compatibility

EMC requirements according to DIN EN 61326-1

Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1: General requirements (IEC 61326-1:2021)

Interference resistance

Suitable for use in industrial areas (Table 2 of the standard)

Interference emission

Class B: Suitable for use in residential areas and areas that are connected to a low voltage network that (also) supplies residential buildings.

NOTE: This equipment has been tested and found to comply with the limits for a Class B digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference in a residential installation. This equipment generates, uses and can radiate radio frequency energy and, if not installed and used in accordance with the instructions, may cause harmful interference to radio communications. However, there is no guarantee that interference will not occur in a particular installation. If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try to correct the interference by one or more of the following measures:

- Reorient or relocate the receiving antenna.
- Increase the separation between the equipment and receiver.
- Connect the equipment into an outlet on a circuit different from that to which the receiver
- Consult the dealer or an experienced radio/TV technician for help.

Shielded Cables: Connections between the devices and peripherals must be made using shielded cables in order to maintain compliance with FCC radio frequency emission limits.

Modifications: Any modifications made to this devices that are not approved by Sartorius may void the authority granted to the user by the FCC to operate this equipment.

14.5 Product Water Quality

14.5.1 Arium® Mini Plus and Arium® Mini Plus | Extend

	Unit	Value (ultrapure water stage)	Value (Pretreatment stage)
Water type		Ultrapure water ASTM Type 1	Pure water type 3
Production output ¹	l/h	–	8
Flow output water dispense ²	l/min	≤ 1.0	Depressurized via ball valve
Volume input ²	l	Between 0.05 and 5 (in 50 ml increments)	–
Typical conductivity	µS/cm	0.055 (compensated at 25 °C ⁵)	< 20 (compensated at 25 °C ⁶)
Typical resistivity	MΩ + cm	18.2 (compensated at 25 °C ⁵)	< 0.05 (compensated at 25 °C ⁶)
TOC content ³ (system with UV lamp)	ppb	≤ 5	–
Content of microorganisms ⁴	CFU/ml	< 0.001	< 0.001
Particle content ⁴	µm	No particles > 0.2 µm	No particles > 0.2 µm
Typical ion retention	%	–	≤ 98
Retention of dissolved organic substances (MW > 300 Dalton)	%	–	> 99
Particle and microorganism retention	%	–	> 99

¹ Depending on the feed water pressure, temperature, and condition of the RO modules

² Depending on the hydrostatic pressure, and connected accessories and final filter

³ Determined with municipal water (Goettingen), TOC approx. 1000 ppb

⁴ When using an Arium® Sterile Plus final filter (Sartopore® 2 150)

⁵ Measured value display at 25 °C compensated or not compensated

⁶ Depending on inlet water

14.5.2 Arium® Mini, Arium® Mini Essential and Arium® Mini Essential | Extend

	Unit	Value
Water type		Ultrapure water ASTM Type 1
Flow output water dispense ²	l/min	≤ 1.0
Volume input ²	l	Between 0.05 and 5 (in 50 ml increments)
Typical conductivity	µS/cm	0.055 (compensated at 25 °C ⁵)
Typical resistivity	MΩ + cm	18.2 (compensated at 25 °C ⁵)
TOC content ³ (system with UV lamp)	ppb	≤ 5
Content of microorganisms ⁴	CFU/ml	< 0.001
Particle content ⁴	µm	No particles > 0.2 µm

² Depending on the hydrostatic pressure, and connected accessories and final filter

³ Determined with municipal water (Goettingen), TOC approx. 1000 ppb

⁴ When using an Arium® Sterile Plus sterile filter (Sartopore® 2 150)

⁵ Measured value display at 25 °C compensated or not compensated

14.6 Feed Water Quality

14.6.1 Arium® Mini Plus and Arium® Mini Plus | Extend

	Unit	Value
Suitability / type		Exclusively potable tap water pursuant to the drinking water standards of the USA, the European Union or Japan.
Inlet pressure	bar	0.5 to 6 (recommended: > 2)
Temperature	°C	2 to 30
TOC	ppb	< 2,000
Max. total hardness (max. CaCO ₃)	ppm	360
Free chlorine	ppm	< 4
Iron (total Fe content)	ppm	< 0.1
Fouling Index (SDI)		< 10
Turbidity	NTU	< 1
pH value		4 to 10

14.6.2 Arium® Mini

	Unit	Value
Suitability / type		Purified water using reverse osmosis, distillation or deionization
Inlet pressure		Depressurized
Temperature	°C	2 to 30
Specific conductivity	µS/cm	<100 (compensated to 25 °C)
TOC	ppb	<50
Turbidity	NTU	<1
pH value		4 to 10

14.6.3 Arium® Mini Essential and Arium® Mini Essential | Extend

	Unit	Value
Suitability / type		Purified water using reverse osmosis, distillation or deionization
Inlet pressure		0 to 6 (recommended: > 2)
Temperature	°C	2 to 30
Specific conductivity	µS/cm	<100 (compensated to 25 °C)
TOC	ppb	<50
Turbidity	NTU	<1
pH value		4 to 10

14.7 Device Properties

	Unit	Value
Devices without a dispensing unit		
Dimensions (width x height x depth)	mm	280 x 510 x 530
Empty weight, approx.	kg	13
Operating weight, approx.	kg	23
Devices with a dispensing unit		
Dimensions (width x height x depth)	mm	370 x 640 x 610
Empty weight, approx.	kg	14
Operating weight, approx.	kg	24
Water treatment method		Adsorption using spherical activated charcoal, catalyst, reverse osmosis, ion exchange, optional UV radiation, and end position particle/sterile filtration or removal of endotoxins, RNases, and DNases.

15 Consumables

This table contains an excerpt of accessories that can be ordered.
For information on other products, contact Sartorius.

Description	Order number
Final filter	
Arium® sterile filter Sterile Plus (Sartopore® 2 150 capsule)	5441307H4--CE
Arium® Cell Plus ultrafilter	H2O-CUF
Arium® pretreatment cartridge ¹	H2O-CPR
Arium® Scientific Pack (ultrapure water cartridge)	H2O-S-PACK
Arium® 5-liter bag ²	H2O-CBS-5-S
Arium® UV lamp	H2O-CEL1

¹ Only required for Arium® Mini Plus

² Only required for Arium® Mini Plus or Arium® Mini

16 Sartorius Service

The Sartorius Service is at your disposal for queries regarding the device.
For information about the service addresses, services provided or to contact a local representative, please visit the Sartorius website (www.sartorius.com).

When contacting Sartorius Service with questions about the system or in the event of malfunctions, be sure to keep the device information—e.g. serial number, hardware, firmware, configuration—close at hand. This information can be found on the manufacturer's ID label and in the "Device Information" menu.

17 Conformity

17.1 EC/EU Declaration of Conformity

The attached Declaration of Conformity hereby confirms compliance of the device with the directives cited.

17.2 UK Declaration of Conformity

The attached Declaration of Conformity hereby confirms compliance of the device with the directives cited.



Original

EG-/EU-Konformitätserklärung EC / EU Declaration of Conformity

Hersteller **Sartorius Lab Instruments GmbH & Co. KG**
 Manufacturer 37070 Goettingen, Germany

erklärt in alleiniger Verantwortung, dass das Betriebsmittel
declares under sole responsibility that the equipment

Geräteart **Reinstwassersystem Arium® Mini Plus, Arium® Mini, Arium® Mini Essential,
 Arium® Mini Plus | Extend, Arium® Mini Essential | Extend**

Device type *Ultrapure water treatment system Arium® Mini Plus, Arium® Mini, Arium® Mini Essential,
 Arium® Mini Plus | Extend, Arium® Mini Essential | Extend*

Baureihe **H2O-MA-T, H2O-MA-T-US, H2O-MA-UV-T, H2O-MA-UV-T-US (Arium® Mini Plus)
 H2O-MM-T, H2O-MM-T-US, H2O-MM-UV-T, H2O-MM-UV-T-US (Arium® Mini)
 H2O-MU-T, H2O-MU-T-US, H2O-MU-UV-T, H2O-MU-UV-T-US (Arium® Mini Essential)
 H2O-MAR-UV-T, H2O-MAR-T (Arium® Mini Plus | Extend)
 H2O-MUR-UV-T, H2O-MUR-T (Arium® Mini Essential | Extend)**

in der von uns in Verkehr gebrachten Ausführung allen einschlägigen Bestimmungen der
 folgenden Europäischen Richtlinien entspricht und die anwendbaren Anforderungen folgender
 harmonisierter Europäischer Normen einschließlich deren zum Zeitpunkt der Erklärung
 geltenden Änderungen erfüllt:

*in the form as delivered fulfils all the relevant provisions of the following European Directives
 and meets the applicable requirements of the harmonized European Standards including any
 amendments valid at the time this declaration was signed listed below:*

Richtlinie
 Directive
 Norm(en)
 Standard(s)

EMV / EMC	RoHS	Maschinen / Machines
2014/30/EU	2011/65/EU	2006/42/EG 2006/42/EC
EN 61326-1:2013 EN IEC 61326-1:2021	EN IEC 63000:2018	EN ISO 12100:2010 EN 61010- 1:2010+A1:2019 ^{*)}

Die Person, die bevollmächtigt ist, die technischen Unterlagen zusammenzustellen:
The person authorized to compile the technical file:

Sartorius Lab Instruments GmbH & Co. KG
 Electronics & Product Compliance
 37070 Goettingen, Germany

Sartorius Lab Instruments GmbH & Co. KG
 Goettingen, 2024-04-30

Mario Hespe
 Managing Director

Halil Yildirim
 Senior Compliance Manager (LPS)

*: angewandte, jedoch für Maschinen nicht harmonisierte Norm / *applied standard, which however is
 not harmonized for machines*

Doc: 2036499-08 SLI15CE018-08.de,en 1 / 1 PMF: 2036498 OP-113_fo1_2020.07.07



Original

SARTORIUS

UK Declaration of Conformity

Manufacturer Sartorius Lab Instruments GmbH & Co. KG
37070 Goettingen, Germany

declares under sole responsibility that the equipment.

Device type Ultrapure water treatment system

Type series Arium® Mini Plus, Arium® Mini, Arium® Mini Essential, Arium® Mini Plus | Extend,
Arium® Mini Essential | Extend

Model H2O-MA-T, H2O-MA-T-US, H2O-MA-UV-T, H2O-MA-UV-T-US (Arium® Mini Plus)
H2O-MM-T, H2O-MM-T-US, H2O-MM-UV-T, H2O-MM-UV-T-US (Arium® Mini)
H2O-MU-T, H2O-MU-T-US, H2O-MU-UV-T, H2O-MU-UV-T-US (Arium® Mini essential)
H2O-MAR-UV-T, H2O-MAR-T (Arium® Mini Plus | Extend)
H2O-MUR-UV-T, H2O-MUR-T (Arium® Mini Essential | Extend)

in the form as delivered fulfils all the relevant provisions of the following British Regulations and meets the applicable requirements of the British Designated Standards including any amendments valid at the time this declaration was signed listed below:

The Electromagnetic Compatibility Regulations 2016

UK Statutory Instruments 2016 No. 1091

BS EN 61326-1:2013; BS EN IEC 61326-1:2021

The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012

UK Statutory Instruments 2012 No. 3032

BS EN IEC 63000:2018

The Supply of Machinery (Safety) Regulations 2008

UK Statutory Instruments 2008 No. 1597

BS EN ISO 12100:2010, BS EN 61010-1:2010+A1:2019*)

The person authorised to compile the technical file:

Sartorius UK Ltd.
Longmead Business Centre, Blenheim Road
KT19 9 QQ Epsom, Surrey, UK

Sartorius Lab Instruments GmbH & Co. KG
Goettingen, 2024-04-30

Mario Hespe
Managing Director

Halil Yildirim
Senior Compliance Manager (LPS)

*: applied standard, which however is not harmonized for machines

Doc: 2726148-02 SLI21UKCA004-02.en 1 / 1 PMF: 2036498 OP-113_fo18_2021.01.01

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany

Phone: +49 551 308 0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote all genders.

Copyright notice:

These instructions, including all components, are protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

Last updated:

04 | 2024