

## Operating Instructions

Original Operating Instructions

# Picus<sup>®</sup> 2

Electronic Pipette

Single-Channel Models | Multi-Channel Models



1000125943



# SARTORIUS



# Contents

<b>1</b>	<b>About This Manual</b> .....	<b>7</b>
1.1	Validity .....	7
1.2	Supporting Documents .....	7
1.3	Target Groups .....	8
1.4	Symbols Used .....	8
1.4.1	Warnings in Operation Descriptions .....	8
1.4.2	Other Symbols Used .....	8
<b>2</b>	<b>Safety Instructions</b> .....	<b>9</b>
2.1	Intended Use .....	9
2.1.1	Modifications to the Device .....	9
2.2	Qualifications of Personnel .....	9
2.3	Functional Capability of the Device Parts .....	9
2.4	Electrical Equipment .....	10
2.4.1	Damage to the Electrical Equipment of the Device .....	10
2.5	Risk of Injury During Continuous Pipetting .....	10
<b>3</b>	<b>Device Description</b> .....	<b>11</b>
3.1	Device Overview .....	11
3.2	Controls and Connection .....	12
3.3	Pipette Tips and Safe-Cone Filters .....	13
3.4	Initial Adjustment and Follow-up Adjustment .....	14
3.5	Sartorius Mobile App .....	14
3.6	Symbols on the Package, on the Device and as Digital Labels .....	14
<b>4</b>	<b>Operating Design</b> .....	<b>16</b>
4.1	Controls .....	16
4.2	Operating Mode Display .....	17
4.2.1	Displays in Menu .....	18
4.3	Messages .....	19
4.3.1	Function of the Softkeys .....	19
4.4	Indicators in the Display .....	20
4.5	Navigating in Menus .....	21
4.5.1	Menus in the Device .....	23
4.6	Menu Structure in the Main Menu .....	23
4.6.1	Menu Structure in the "Settings" Sub-menu .....	24
4.7	Menu Structure in the "Advanced Functions" Menu .....	25
4.8	Menu Structure in the "Memory Places" Menu .....	25
4.9	List of Parameters .....	26
4.9.1	Parameters in the "Advanced Functions" Menu .....	26
4.9.2	Parameters in the "Settings"/"Languages" Menu .....	27
4.9.3	Parameters in the "Settings"/"Theme" Menu .....	27
4.9.4	Parameters in the "Settings"/"Bluetooth" Menu .....	27
4.9.5	Parameters in the "Settings"/"Password" Menu .....	27
4.9.6	Parameters in the "Settings"/"Sound" Menu .....	28
4.9.7	Parameters in the "Settings"/"Backlight" Menu .....	28
4.9.8	Parameters in the "Settings"/"Date & time" Menu .....	28
4.9.9	Parameters in the "Settings"/"Tip Ejection" Menu .....	28

4.9.10	Parameters in the “Settings”/“Adjustment” Menu	29
4.9.11	Parameters in the “Settings”/“User ID” Menu	29
4.9.12	Parameters in the “Settings”/“Reminder” Menu	30
4.9.13	“Settings”/“Reset” Menu	30
<b>5</b>	<b>Installation</b>	<b>31</b>
5.1	Scope of Delivery	31
5.2	Unpacking	31
5.3	Acclimatization	31
5.4	Inserting or Removing a Safe-Cone Filter	32
5.4.1	Inserting a Safe-Cone Filter	32
5.4.2	Removing the Safe-Cone Filter	32
<b>6</b>	<b>Commissioning</b>	<b>33</b>
6.1	Clean the Device	33
6.2	Charging the Device	33
6.2.1	Charging with the Charging Carousel or Charging Stand	33
6.2.2	Charging with the USB Charger	33
6.3	Switching the Device On or Off	34
6.3.1	Returning to Active Mode	34
6.4	Rinsing Tips	34
<b>7</b>	<b>System Settings</b>	<b>35</b>
7.1	Setting Defaults	35
7.2	Using the Device with the Sartorius Pipetting MobileApp	35
7.2.1	Downloading the Sartorius Pipetting MobileApp	35
7.2.2	Adding the Device in the app	35
7.2.3	Creating the Pipette Set	36
7.2.4	Starting the Workflow	36
7.3	Performing a Software Update	36
7.4	Managing User Data	37
7.4.1	Creating a User ID	37
7.4.2	Activating Password Protection	37
7.4.3	Logging Into or Out of the Device	38
7.4.4	Deactivating Password Protection	38
7.5	Adjusting the Device	39
7.5.1	Activating the Adjustment Setting	41
7.5.2	Documenting Adjustment for Liquids Other Than Water	41
7.6	Setting Up Reminders	41
7.6.1	Deactivating Reminders	42
<b>8</b>	<b>Operation</b>	<b>43</b>
8.1	Accessing Digital Compliance Labels	43
8.2	Using Operating Modes (Examples)	43
8.2.1	Multi Dispensing	43
8.2.2	Manual Pipetting	44
8.3	Locking and Unlocking the Device	45
8.3.1	Activating the Device Lock	45
8.3.2	Opening the Device Lock	45
<b>9</b>	<b>Cleaning and Maintenance</b>	<b>46</b>
9.1	Cleaning the Surfaces of the Device	46
9.2	Maintenance Schedule	46

9.3	Cleaning and Greasing the Lower Assembly of the Device	47
9.3.1	Preparing the Device	47
9.3.2	Dismantling the Lower Assembly (1-channel Models)	47
9.3.3	Cleaning Individual Components (1-channel Models)	49
9.3.4	Greasing Components (1-channel models)	50
9.3.5	Assembling the Lower Assembly (1-channel Models)	51
9.4	Sterilising the Device	52
9.4.1	Selecting the Sterilization Method	52
9.4.2	Sterilizing the Device with UV Radiation	52
9.4.3	Sterilizing the Device with an Decontamination Agent	52
9.4.4	Sterilising the Lower Assembly of the Device	53
9.5	Defining the Test Routine for the Routine Test	54
9.6	Performing the Routine Test	56
9.6.1	Aspirating the Test Water	56
9.6.2	Dispensing Test Water and Performing the Measurement	57
9.6.3	Evaluating Measured Values	57
<b>10</b>	<b>Trouble</b>	<b>59</b>
10.1	Troubleshooting	59
10.2	Resetting to Defaults	60
<b>11</b>	<b>Decommissioning</b>	<b>61</b>
11.1	Decommissioning the Device	61
<b>12</b>	<b>Storage and Shipping</b>	<b>61</b>
12.1	Storage	61
12.2	Returning the Device and Parts	62
<b>13</b>	<b>Disposal</b>	<b>62</b>
13.1	Disposing of the Device and Parts	62
<b>14</b>	<b>Technical Data</b>	<b>63</b>
14.1	Dimensions and Weights	63
14.1.1	1-channel models	63
14.1.2	Multi-channel models with 8	63
14.1.3	Multi-channel models with 12	63
14.2	Ambient Conditions	64
14.3	Electrical Data	64
14.3.1	Power Supply	64
14.3.2	Safety of Electrical Equipment	65
14.3.3	Safety and Electromagnetic Compatibility	65
14.4	Interfaces	65
14.5	Operating Modes with Advanced Functions	66
14.6	Acceleration Sensor	66
14.7	RFID Tag	66
14.8	Adjustment	67
14.8.1	Initial Adjustment	67
14.9	Materials	67
14.10	Cleaning and Sterilization	68
14.10.1	Cleaning Agents, Decontamination Agents and Cleaning Methods	68
14.10.2	Approved Sterilization Methods for Lower Assembly	68
14.11	Battery	69

14.12	Routine test.....	69
14.12.1	Conditions for Aspirating the Test Liquid.....	69
14.12.2	Correction Factor Z ( $\mu\text{L} / \text{mg}$ ) for Routine Test.....	69
14.13	Performance Specification.....	70
14.13.1	Single-Channel Models.....	70
14.13.2	Multi-Channel Models.....	72
14.14	Speed Table.....	74
14.14.1	1-channel Pipette.....	74
14.14.2	Multi-channel Pipette.....	74
<b>15</b>	<b>Accessories and Consumables.....</b>	<b>75</b>
15.1	Accessories.....	75
15.2	Consumables.....	75
<b>16</b>	<b>Sartorius Service.....</b>	<b>76</b>
<b>17</b>	<b>Conformity of the Device.....</b>	<b>76</b>
17.1	Compliance Information and Information for Users for Radio Equipment.....	78
<b>18</b>	<b>Trademark Information.....</b>	<b>78</b>
18.1	Bluetooth®.....	78

# 1 About This Manual

## 1.1 Validity

This manual is part of the device and must be read in full and kept safe. The manual applies to the following versions of the device:

Device	Volume range ( $\mu\text{L}$ )	Maximum volume / nominal volume ( $\mu\text{L}$ )	Item number
Picus <sup>®</sup> 2 electronic pipette			
1-channel model	0.5 - 10	10	LH-747021
	5 - 120	120	LH-747041
	10 - 300	300	LH-747061
	50 - 1000	1000	LH-747081
	100 - 5000	5000	LH-747101
	500 - 10000	10000	LH-747111
8-channel model	0.5 - 10	10	LH-747321
	5 - 120	120	LH-747341
	10 - 300	300	LH-747361
	50 - 1200	1200	LH-747391
12-channel model	0.5 - 10	10	LH-747421
	5 - 120	120	LH-747441
	10 - 300	300	LH-747461
	50 - 1200	1200	LH-747491

## 1.2 Supporting Documents

- ▶ In addition to this manual, please note the following documentation:
  - Instructions for the consumables, accessories, e.g. pipette stands, pipette tips

## 1.3 Target Groups

This manual is aimed at the following target groups. The target groups must have the specified knowledge.

Target group	Knowledge and qualifications
Operator	The operator is familiar with the device and the associated work processes. The operator is aware of the dangers that can occur when working with the device, and can avoid these dangers.*
Administrator	The administrator is responsible for integrating the device into the production process. The administrator ensures reliable operation of the system and device software.*

\* When a person from the target group operates the software interface of the device, they are simultaneously a "user".

## 1.4 Symbols Used

### 1.4.1 Warnings in Operation Descriptions

#### CAUTION

Identifies a hazard that can result in moderate or minor injuries if **not** avoided.

### 1.4.2 Other Symbols Used

- ▶ Action: describes activities that must be executed. The activities in action sequences must be executed in consecutive order.
- ▷ Result: describes the result of the activities executed.
- [ ] Refers to controls and display elements. Identifies status messages, warning messages and error messages.

#### Illustrations in this manual

Depending on the device configuration, the illustrations of the device and user display may differ slightly from the supplied device. The variants shown in this manual are examples.



## 2 Safety Instructions

### 2.1 Intended Use

The Picus® 2 pipette is a general purpose laboratory device. The device is intended, designed and manufactured for dispensing liquids in a variety of applications.

The device's volume range must be taken into consideration when selecting the application (see Chapter "14.13 Performance Specification", page 70). Volume deviations can be caused by dosing liquids with different physical properties.

Pipette tips (tips) must be used with the device. Sartorius recommends using only Sartorius tips (see Chapter "15.2 Consumables", page 75).

The tips are designed for single use and must be disposed of after being used once.

The device is exclusively intended for use in accordance with these instructions. Any other use is considered **unintended** use.

#### Usage conditions for the device

Do **not** use the device in potentially explosive environments. Only use the device indoors.

Only use the device with the equipment and under the operating conditions described in the technical data in this manual.

#### 2.1.1 Modifications to the Device

If the device is modified: People could be put at risk. Device-specific documents and product approvals could lose their validity.

If you have any queries regarding modifications to the device, please contact Sartorius.

### 2.2 Qualifications of Personnel

People without adequate knowledge of how to handle the device safely could injure themselves and others.

If a specific qualification is required for an activity: The target group is specified. If **no** qualification is specified: The activity can be executed by the "Operator" target group.

### 2.3 Functional Capability of the Device Parts

Device parts that are **not** functional, e.g. due to damage or wear, can cause malfunctions. This could cause injury to people.

► If device parts are **not** functional: Do **not** use the device.

## 2.4 Electrical Equipment

### 2.4.1 Damage to the Electrical Equipment of the Device

Damage to the electrical equipment of the device, e.g. damage to the insulation, can lead to unforeseeable hazards.

- ▶ If there are defects in the electrical equipment, disconnect the device from the power supply and contact Sartorius Service.
- ▶ Keep moisture away from live parts. The moisture can cause short circuits.

## 2.5 Risk of Injury During Continuous Pipetting

An unsuitable posture during pipetting, or using the device without taking breaks, can result in musculoskeletal disorders or repetitive strain injury (RSI) to the hand.

- ▶ When pipetting, take regular breaks and relax your hand.
- ▶ When pipetting, stand or sit upright.
- ▶ Place your index finger under the hook and operate the device with your thumb.
- ▶ Do **not** tense up while pipetting.

## 3 Device Description

### 3.1 Device Overview

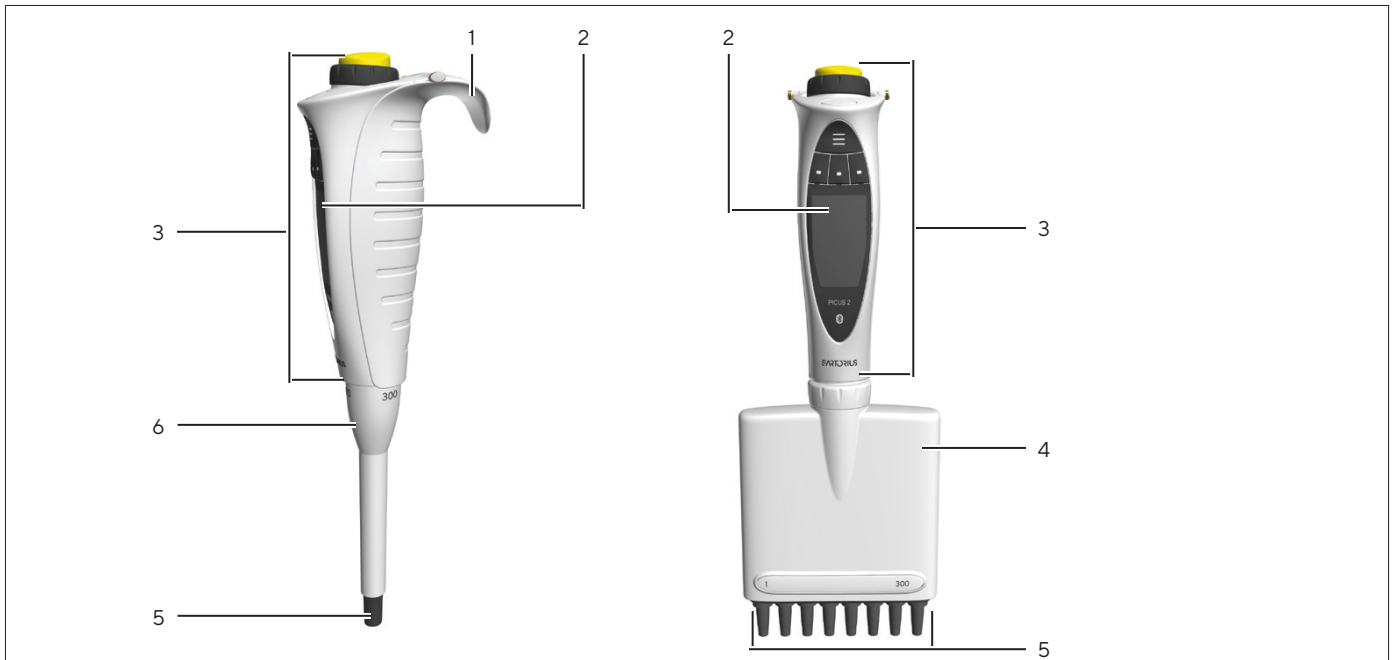


Fig. 1: 1-channel model and multi-channel model (example)

Pos.	Name	Description
1	Finger Hook	For resting the device on your finger.
2	User display	Displays all the relevant information.
3	Control head	To control the device and trigger individual pipetting steps
4	Dispensing head	Consists of: <ul style="list-style-type: none"> <li>– Piston</li> <li>– Cylinder</li> <li>– Tip cone</li> <li>– Tip ejection system</li> </ul>
5	Tip Cone(s)	<ul style="list-style-type: none"> <li>– For attaching tips</li> <li>– Optiload system to load tips evenly (only for multi-channel models)</li> </ul>
6	Ejection sleeve	<ul style="list-style-type: none"> <li>– Electronically ejects the tip(s).</li> <li>– With 1-channel models: Height-adjustable</li> <li>– Contains the specification of the maximum volume.</li> </ul>

## 3.2 Controls and Connection

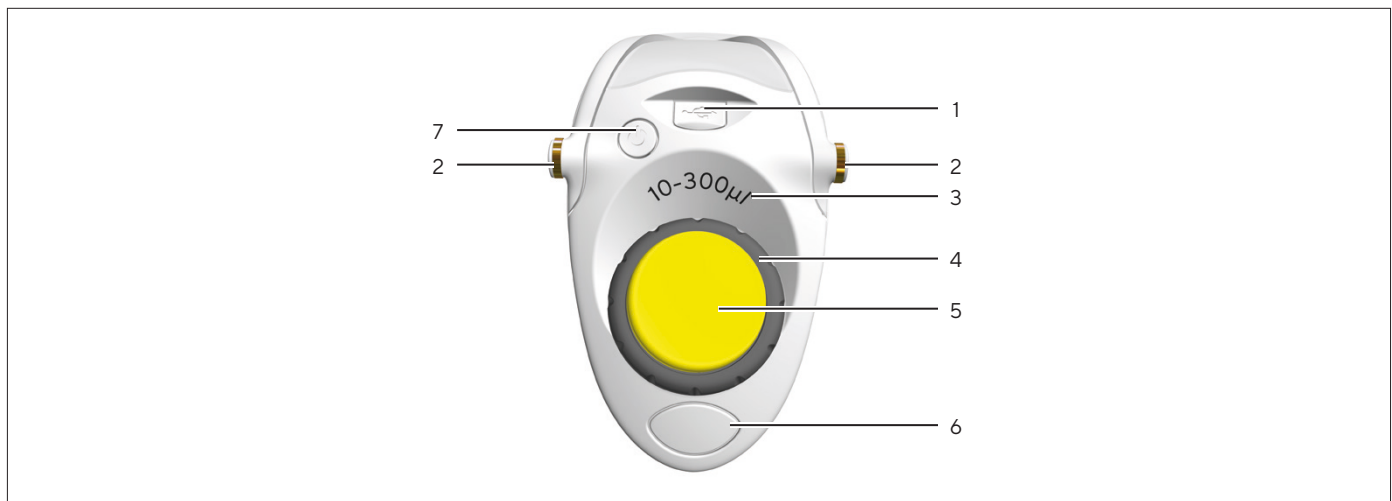


Fig.2: Controls and connection (example)

Pos.	Name	Description
1	Micro USB port	For connection to the mains plug or a to PC
2	Charging contact	Is intended to charge the device with the charging stand or carousel.
3	Volume range	Indicates the volume range.
4	Scrollwheel	<ul style="list-style-type: none"> <li>– Allows you to scroll through menus and set values.</li> <li>– Controls the aspiration and dispensing of liquids in manual pipetting mode.</li> </ul>
5	Operating button	<ul style="list-style-type: none"> <li>– Used to confirm settings and trigger the piston movement during aspiration, dispensing and repeated blow-out.</li> <li>– The colour indicates the volume range (for the colour of the operating button, see Chapter “14.13 Performance Specification”, page 70).</li> </ul>
6	Electronic tip ejection	Is intended to eject the tip.
7	On   Off button	For switching the device on or off

### 3.3 Pipette Tips and Safe-Cone Filters

The pipette tip must be adapted to the volume range of the device and the degree of purity. The colour of the operating button and the colour of the tip trays must match. We recommend using Sartorius Optifit Tips or Safetyspace® Filter Tips. The use of Safe-Cone Filters or Safetyspace® Filter Tips helps to prevent contamination of the device. Safe-cone filters are replacable and can be used together with Optifit tips. Safe-Cone filters are inserted in pipette tip cones.

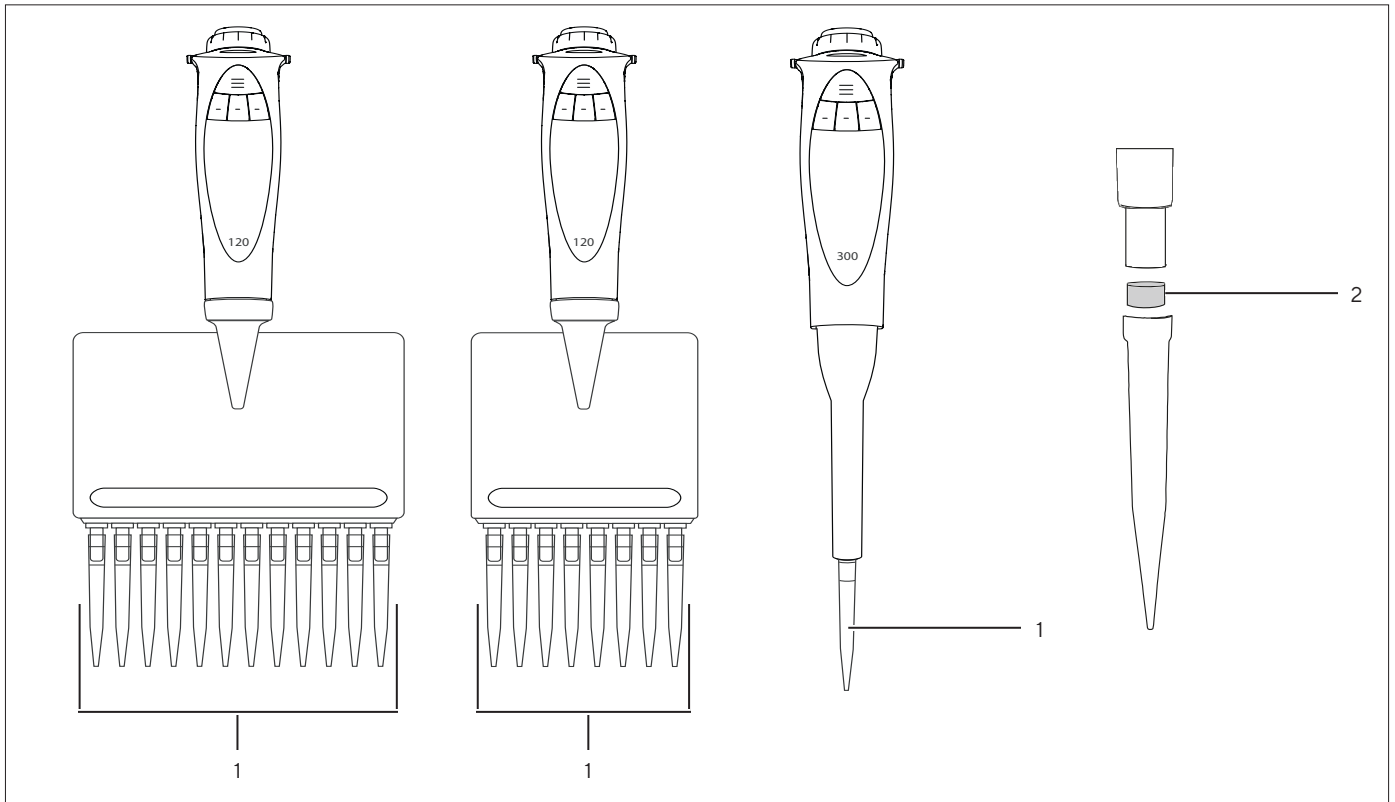


Fig. 3: Pipette tips on multi-channel models and 1-channel model and Safe-Cone-Filter (example)

Pos.	Name	Description
1	Pipette tips	Serve to aspirate and dispense liquids.
2	Safe-Cone-Filter	<ul style="list-style-type: none"> <li>– Helps prevent liquids from contaminating the device.</li> <li>– Available separately for models over 10 µl</li> </ul>

## 3.4 Initial Adjustment and Follow-up Adjustment

Initial adjustment: The device is adjusted on delivery in accordance with ISO 8655-1 (see Chapter “14.8.1 Initial Adjustment”, page 67).






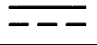
Follow-up adjustment: The device can be re-adjusted, e.g., if it is determined that the delivered volume is **not** within the maximum number of permissible errors or a liquid other than water is to be pipetted. The described procedure must be used for the adjustment. The current adjustment of the device is displayed in the “Settings / Adjustment” menu.

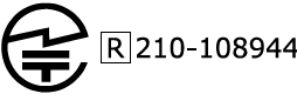







## 3.5 Sartorius Mobile App

The device can also be used with companion app (Chapter “7.2 Using the Device with the Sartorius Pipetting MobileApp”, page 35). In the app you will have multiple options to speed up your routine workflows and customize your pipettes. The app can also be used to perform device updates (Chapter “7.2.3 Creating the Pipette Set”, page 36). The app is available at App store and Play store.



## 3.6 Symbols on the Package, on the Device and as Digital Labels

Symbol	Meaning
	Bluetooth® symbol: Indicates that a Bluetooth connection can be established with this device.
	Autoclaving symbol: Component is autoclavable.
	The device complies with applicable EU directives and standard.
	The device complies with applicable UK legislation and standards.
	Separate collection of electronic equipment
	Direct current

Symbol	Meaning
	GITEKI symbol: The device complies with the standards and regulations of the Japanese Radio Law.
	The device complies with the Official Mexican Standards.
	This device contains certain hazardous substances and can be used safely during its environmental protection use period.
	The device meet, along with the approved stipulated frequency, the strict Radio Frequency Interference standards stipulated by South Africa's independent communications authority (ICASA).
	The device complies with the legal requirements for radio technology products in the Taiwanese market.
	The device complies with the Korean Radio Waves Act.
<b>R-NZ</b>	The device complies with applicable New Zealand radio product rules.
	The device complies with the U.S. Federal Communications Commission (FCC) regulations.
	The device complies with applicable Australian rules.

## 4 Operating Design

### 4.1 Controls

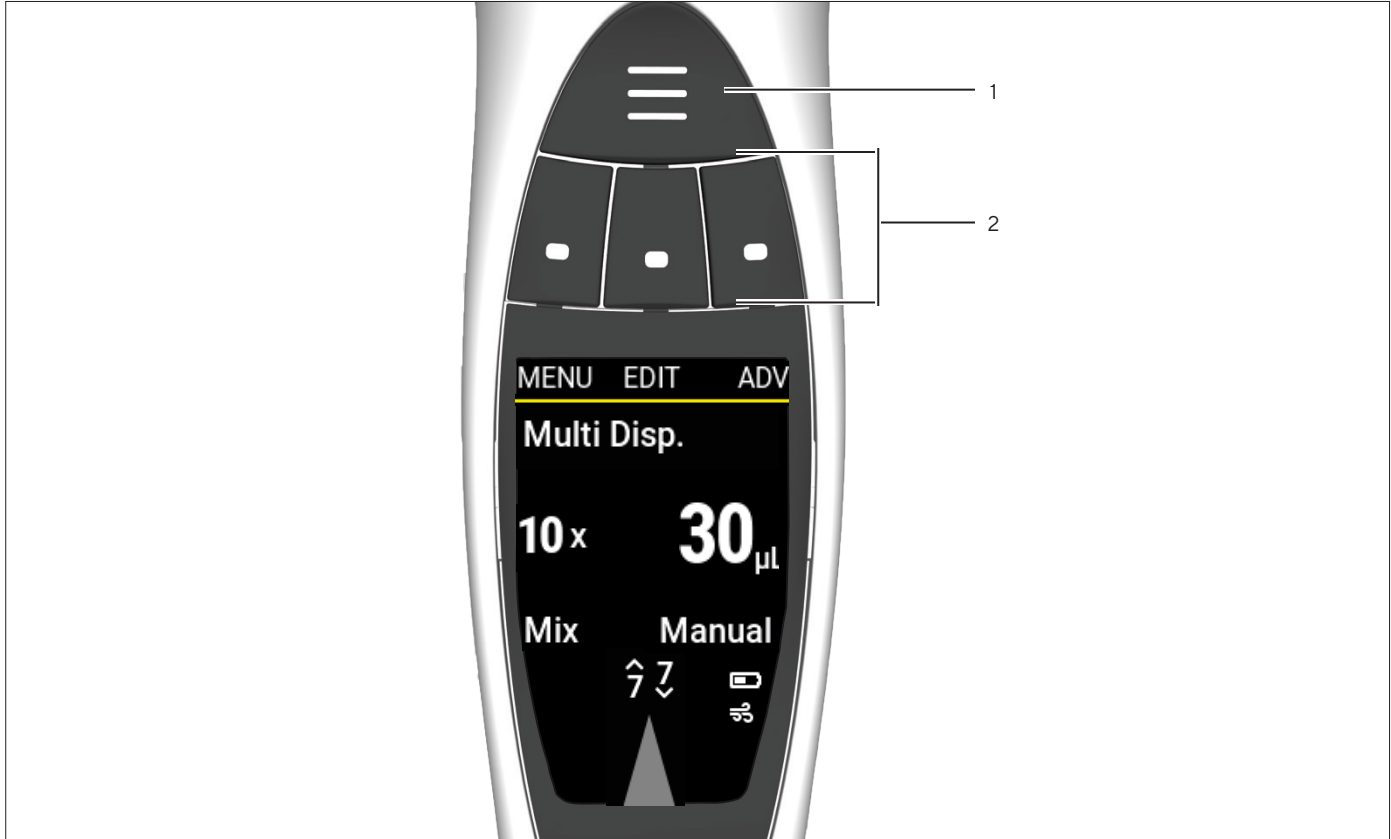


Fig.4: Controls on the device

Pos.	Name	Description
1	Hotkey	Save and activate frequently used or favorite pipette settings.
2	Softkeys	<ul style="list-style-type: none"> <li>– For controlling the device.</li> <li>– The functions of the individual keys vary according to the display.</li> </ul>



## 4.2 Operating Mode Display

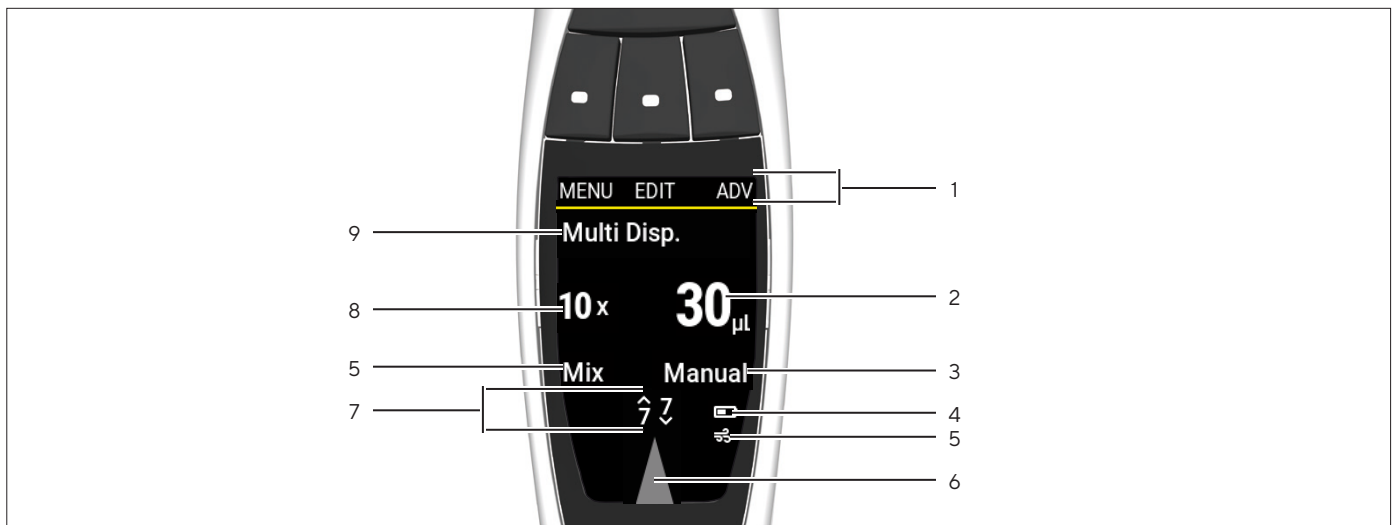


Fig. 5: Operating mode display (example)

Pos.	Name	Description
1	Softkey function	Indicates which function is assigned to the overlying softkey.
2	Pipetting volume	<ul style="list-style-type: none"> <li>– Indicates what volumes are being taken up or dispensed..</li> <li>– In Multi Dispense mode: Indicates partial volumes dispensed per dispensing process.</li> </ul>
3	Advanced functions property	Adjustable parameters for the advanced functions. Value and unit vary according to the activated advanced function.
4	Battery indicator	Displays the battery level.
5	Advanced functions indicator	Displays the activated advanced function.
6	Aspirate   dispense display arrow	<ul style="list-style-type: none"> <li>– If the tip is pointing up: The device is set up to aspirate</li> <li>– If the tip is pointing down: The device is set up to dispense.</li> </ul>
7	Speed	Indicates the speed with which the device aspirates or dispenses.
8	Pipetting activity number	Indicates how often aspiration   dispensing is carried out in the current operating mode. This parameter is only available in the multi-dispensing, sequentiell dispensing and multi-aspiration operating modes.
9	Operating mode	Indicates the selected operating mode.

### 4.2.1 Displays in Menus

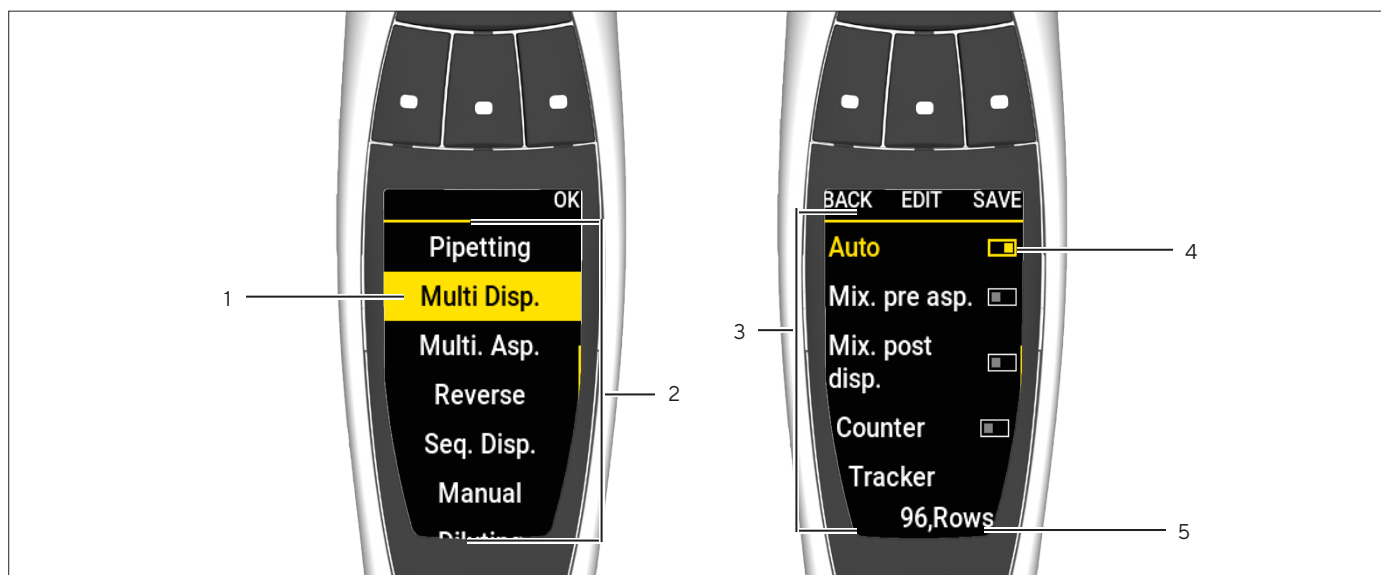


Fig. 6: Displays in menus (example)

Pos.	Name	Description
1	Menu entry	Yellow background: Currently selected.
2	Main menu	Contains available operating modes and settings menu.
3	Advanced functions	Number and type of advanced functions. The number varies according to the activated operating mode.
4	On   Off bar	<ul style="list-style-type: none"> <li>– If the dark bar is on the right: The function is activated.</li> <li>– If the dark bar is on the left: The function is deactivated.</li> </ul>
5	Selection parameters	Are Predefined and selectable with the scrollwheel.

## 4.3 Messages

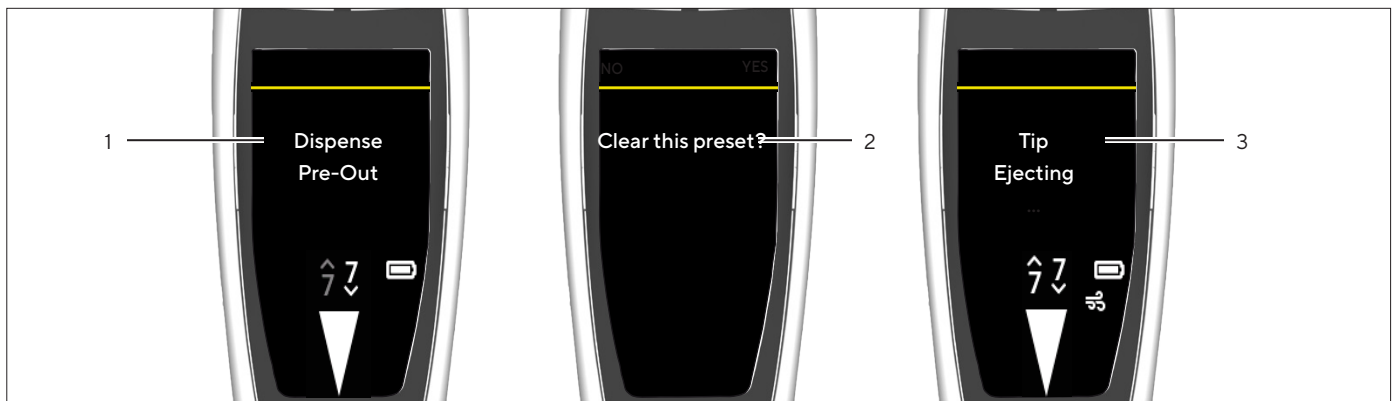











Fig.7: Messages (example)











Pos.	Name	Description
1	Prompt	<ul style="list-style-type: none"> <li>– Shows what the device will do next.</li> <li>– Initiation via the operating button is required.</li> </ul>
2	Query	<ul style="list-style-type: none"> <li>– Shows the operator options.</li> <li>– The selection is made using the softkeys.</li> </ul>
3	Notification	<ul style="list-style-type: none"> <li>– Shows what the device is doing.</li> <li>– <b>No</b> action by the operator is required.</li> </ul>

### 4.3.1 Function of the Softkeys

Symbol	Designation	Description
<b>MENU</b>	[MENU] softkey	Opens the main menu.
<b>ADV</b>	[ADV] softkey	Opens the window with the available advanced functions.
<b>EDIT</b>	[EDIT] softkey	Activates edit mode.
<b>OK</b>	[OK] softkey	<ul style="list-style-type: none"> <li>– In the menu: Opens the selected menu entry.</li> <li>– In edit mode: Applies the selection.</li> <li>– For advanced functions: <ul style="list-style-type: none"> <li>– Opens parameter entries</li> <li>– Activates edit mode</li> <li>– Returns to the menu</li> </ul> </li> </ul>
<b>BACK</b>	[BACK] softkey	Returns to the previous display.
<b>Pick</b>	[Pick] softkey	Selects the current entry.
<b>NEXT</b>	[NEXT] softkey	Skips to the next parameter.
<b>END</b>	[END] softkey	Cancel the pipetting process.
<b>SAVE</b>	[SAVE] softkey	Saves the current program at the selected memory place.
<b>RESET</b>	[RESET] softkey	Resets the date and cycle counter.

Symbol	Designation	Description
	[NO] softkey	<ul style="list-style-type: none"> <li>– Does <b>not</b> apply the changes.</li> <li>– Does <b>not</b> execute the action.</li> </ul>
	[YES] softkey	<ul style="list-style-type: none"> <li>– Applies the changes.</li> <li>– Does execute the action.</li> </ul>
	[PREV] softkey	With activated protocol: Returns to the previous protocol step.
	[NEXT] softkey	With activated protocol: Skips to the next protocol step.
	[ABC] softkey	<ul style="list-style-type: none"> <li>– Indicates that upper case is activated.</li> <li>– With softkey pressed: Switches to [abc] function.</li> </ul>
	[abc] softkey	<ul style="list-style-type: none"> <li>– Indicates that lower case is activated.</li> <li>– With softkey pressed: Switches to [123] function.</li> </ul>
	[123] softkey	<ul style="list-style-type: none"> <li>– Indicates that numbers are activated.</li> <li>– With softkey pressed: Switches to [#@!] function.</li> </ul>
	[#@!] softkey	<ul style="list-style-type: none"> <li>– Indicates that special characters are activated.</li> <li>– With softkey pressed: Switches to [CLEAR] function.</li> </ul>
	[CLEAR] softkey	<ul style="list-style-type: none"> <li>– Deletes the text.</li> <li>– With softkey pressed: Switches to [ABC] function</li> </ul>

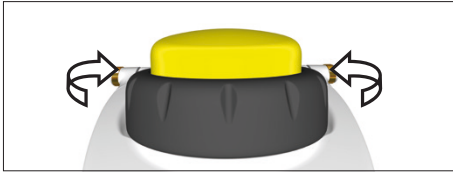
## 4.4 Indicators in the Display

Symbol	Designation	Description
	Indicator [Charge status]	<ul style="list-style-type: none"> <li>– Shows the current charge status of the device.</li> <li>– If the battery level is low: The indicator flashes.</li> </ul>
	Indicator [Indicator arrow]	Indicates which pipetting process is selected: Aspirate or dispense.
	Indicator [Dispensing speed]	Shows the dispensing speed.
	Indicator [Aspiration speed]	Shows the aspiration speed.
	Indicator [Blowout]	Shows that the additional function [Blowout] is activated.
	Indicator [Bluetooth]	Shows that Bluetooth is activated.
	Indicator [Counter]	Shows that the additional function [Counter] is activated.
	Indicator [Plate Tracker]	Shows that the additional function [Plate Tracker] is activated.
	Indicator [Mix]	Shows that the additional function [Mix. pre. asp.] and/or [Mix. post disp] is activated.
	Indicator [Delay]	Shows that the additional function [Auto] is activated.

## 4.5 Navigating in Menus

### Procedure

#### MENU



- ▶ To call up the main menu: Press the [MENU] softkey.
- ▶ When scrolling in menus: Turn the scrollwheel in the desired direction.
  - Counterclockwise: Scroll down.
  - Clockwise: Scroll up.

- ▶ To select a menu option or save settings:
  - ▶ Press the operating button (1) or the [OK] softkey.

- ▶ To activate edit mode for the pipetting settings: Use one of the following options:

- ▶ Turn the scrollwheel as far as it will go in one direction.
- ▶ Press the [EDIT] softkey.

- ▷ Edit mode is activated.

- ▶ To change a parameter: Turn the scrollwheel in the desired direction:
  - Counterclockwise: Increase the value or activate the functions.
  - Turn clockwise: Reduce the value or deactivate the functions.

#### NEXT

- ▶ To apply the changes in edit mode and switch to the next parameter: press the [NEXT] softkey.
- ▷ The next parameter is activated for editing. If there are **no** more parameters, the cursor returns to the first parameter.

#### BACK

- ▶ To return to the next highest menu level or to exit applications without saving: Press the [BACK] softkey.
- ▶ To trigger the piston movement for aspiration, dispensing and repeated blow-out: Press the operating button.
- ▶ To move the piston in manual and titration mode for aspiration and titration: Turn the scrollwheel.

#### ABORT

- ▶ To cancel a mixing operation prematurely: Press the [ABORT] softkey.

#### ADV

- ▶ To call up the advanced functions menu (only possible from a operating mode): Press the [ADV] softkey.



#### OK

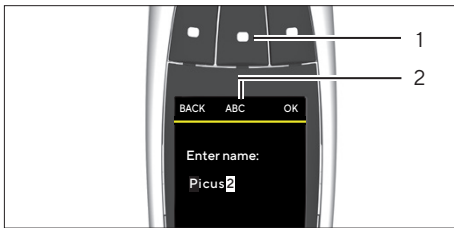
- ▶ To activate a saved program: Press the [OK] softkey.

#### SAVE

- ▶ To save the current program At the selected memory slot press the [SAVE] softkey.

## Text input

- ▶ When entering text, navigate between the positions, e.g. to create the user ID:
  - ▶ In view mode: For one position to the left, turn the scrollwheel anticlockwise, and for one position to the right, turn the scrollwheel clockwise.
  - ▶ In edit mode: Press the [OK] softkey.
  - ▷ The cursor moves one position further to the left.
- ▶ To change an entry during text input:
  - ▶ Activate edit mode.
  - ▶ Turn the scrollwheel until the desired letter appears.
  - ▶ Press the [OK] softkey or the operating button.
  - ▷ The selected letter is applied and the cursor moves one place to the left.
  - ▶ Repeat the settings in this way until the cursor has reached the last position.
  - ▶ To apply the entry: Press the [OK] softkey.
  - ▷ The changed entry is saved and the control returns to the next higher menu level.



- ▶ To enter lower case letters, numbers or symbols during text input: Press the middle softkey (1) until the desired row of characters (2) appears.
- ▶ To delete a position during text input:
  - ▶ Press the middle softkey (1) until the [CLEAR] softkey function appears.
  - ▶ Press the [OK] softkey.
  - ▷ The entry at this position is deleted and the cursor moves one position to the right.

## 4.5.1 Menus in the Device

► Navigating in menus (see Chapter “4.5 Navigating in Menus”, page 21).

Level 1	Level 2	Description
Device	Main menu	<ul style="list-style-type: none"> <li>– Select the operating mode.</li> <li>– Make settings on the device.</li> </ul>
	Advanced functions	Activate advanced functions for a operating mode.
	Memory slots	Save and activate recently used or preferred pipetting settings.

## 4.6 Menu Structure in the Main Menu

Level 1	Level 2	Description
Menu	Pipetting	<p>Aspirate a selected volume of liquid into the tip and then dispense it. This mode is suitable for the following liquids:</p> <ul style="list-style-type: none"> <li>– Aqueous liquids</li> <li>– Liquids with a low quantity of detergents or proteins</li> <li>– Solvents</li> </ul>
	Multi-Disp. “Multi-Dispensing”	<p>Aspirate the total volume and an excess volume of liquid and dispense repeatedly in equal sub-volumes. This mode is suitable for long pipetting rows and for dispensing into microplates.</p>
	Multi. Asp. “Multi-Aspiration”	<p>Set the liquid volume and number of aspiration processes in advance. Then aspirate the selected liquid volumes several times until the cycle is complete. Finally, dispense the entire aspirated liquid volumes in one go. This mode is suitable for merging samples and washing microplates.</p>
	Reverse	<p>Aspirate a selected and an excess volume of liquid. The excess remains in the tip and is disposed of when the tip is ejected. This mode is suitable for the following liquids:</p> <ul style="list-style-type: none"> <li>– Biological liquids</li> <li>– Foaming liquids</li> <li>– Viscous liquids</li> </ul>
	Seq. Disp. “Sequential Dispensing”	<p>Repeated dispensing of selected volumes of liquid in a desired sequence. The liquid volumes can be set directly in the operating mode (1 - 9). This mode is suitable for dilution rows or for creating calibration curves.</p>
	Manual	<p>When aspirating and dispensing liquid volumes, control the piston movement manually with the scrollwheel. This mode is suitable for the following applications:</p> <ul style="list-style-type: none"> <li>– Measuring reagents</li> <li>– Applications in which the pipetting speed must be controlled manually.</li> </ul>
	Diluting	<p>Aspirate the volumes of liquid and dispense at the same time. First the dilution solution is aspirated, then an air cushion and finally the sample or reagent, to avoid contamination. This mode is ideal for e.g. diluting samples and reagents.</p>

Level 1	Level 2	Description
Menu	Titration	Aspirate the entire liquid volumes and determine the dispensing speed manually. The user display shows the dispensed volume in real-time during dispensing.  This mode is suitable for determining the unknown concentration of an identified analyte.
	Settings	Make settings on the device.
	Compliance	Access information on compliance with country-specific directives and regulations and the associated digital labels, e.g. ANATEL or FCC.

#### 4.6.1 Menu Structure in the “Settings” Sub-menu

Level 1	Level 2	Description
Settings	Languages	Set the menu language of the display.
	Themes	Set the UI color by selecting a theme. Default theme selection is model-specific, matching the color of the model’s operating button.
	Bluetooth	Manage the Bluetooth connection.
	Password	Activation of password protection and input of the password. Passwords can be created for an administrator and for a user.
	Sound	Set the sounds on the device.
	Backlight	Set the display lighting.
	Date & time	Enter the date and time.
	Tip ejection	Set the tip ejection settings.
	User ID	Define the user ID. This is displayed on the display when the device is switched on.
	Adjustment	Activate customer-specific adjustment and set the device to the adjustment points.
	Reminder	Set the reminder function for calibration, maintenance, and quick check.
	Information	View the model variant, the current software and Bluetooth® version, and the battery charge level.
	Reset	Reset all modified settings and saved programs to the defaults.
Settings	Pip.lock	Lock the device until the next maintenance measure or cleaning process. The lock prevents any use of the device and the message “Pipette locked, maintenance required” is displayed. We recommend this function when there are device errors or the device is contaminated.



## 4.7 Menu Structure in the “Advanced Functions” Menu

Level 1	Level 2	Description
Advanced functions*	Counter	The counter counts the pipetting cycles.
	Mix. pre. asp.	Activates mixing before aspiration.
	Mix. post disp.	Activates mixing post dispense.
	Auto pre-out	When “Auto pre-out” is activated: The device automatically dispenses pre-out volume before the actual dispensing process is started.
	Excess Adj.	<ul style="list-style-type: none"> <li>– Is intended to define the excess volume (safety volume).</li> <li>– If no excess volume is set: The default value is used.</li> </ul>
	Plate Tracker	<ul style="list-style-type: none"> <li>– The tracker makes it easier to dispense into the correct wells of a microplate by specifying the next position on a microplate during dispensing.</li> <li>– Works with 96- and 384 well plates.</li> </ul>
	Auto	<ul style="list-style-type: none"> <li>– Auto dispensing is intended for automatic dispensing of partial volumes without having to press the operating button.</li> <li>– The dispensing interval must be defined.</li> </ul>
	Fast Vol.	<ul style="list-style-type: none"> <li>– Works with the “Titration” operating mode.</li> <li>– With fast dispensing, a selected quantity of the total volume is dispensed automatically. The remaining volume is then dispensed manually.</li> </ul>
	Blowout	<ul style="list-style-type: none"> <li>– Activates repeated blow-out.</li> <li>– This is used for liquid residues in the filter tip or when dispensing liquids that can leave residues in the filter tip.</li> <li>– Can be combined with other advanced functions.</li> </ul>

\* The “Plate Tracker” and “Counter” advanced functions **cannot** be activated at the same time. The [Blow-out] advanced function can be activated at the same time as all other advanced functions.

## 4.8 Menu Structure in the “Memory Places” Menu

Level 1	Level 2	Description
Memory slots	M1-M20	Is intended to save and activate pipetting settings. Can be accessed via hotkey.

## 4.9 List of Parameters

### 4.9.1 Parameters in the “Advanced Functions” Menu

Parameter	Settings	Settings	Explanation
Counter	ON	0-999	Activates “Counter”. Counter wraps back to 0 if 999 is exceeded.
	OFF*		Deactivates the “Counter” advanced function.
Mix. pre.asp.	ON	Manual Cyclic (1 - 99)	Activates the pre-aspiration mixing function. When activated, it is possible to choose between manual and cyclic mixing.
	OFF*		Deactivates the pre-aspiration mixing function.
Mix. post disp.	ON	Manual Cyclic (1 - 99)	Activates the post-dispensing mixing function. When activated, it is possible to choose between manual and cyclic mixing.
	OFF*		Deactivates the post-dispensing mixing function
Auto pre-out	ON		Activates the automatic dispensing of pre-out volume.
	OFF*		Deactivates the automatic dispensing of pre-out volume.
Excess Adj.	1-25		Enter the safety volume.
Tracker	ON	96, rows 96, columns 384, rows 384, columns	<ul style="list-style-type: none"> <li>– Activates the “Tracker” advanced function.</li> <li>– Set the number of wells.</li> <li>– Set the direction (pipetting vessel in rows   pipetting in columns).</li> <li>– Set the first dispensing position (A1-H12).</li> </ul>
	OFF*		Deactivates the “Tracker”.
Auto	ON	0.0 s-9.9 s	Activates “Auto”. Set the interval until the next dispensing step.
	OFF*		Deactivates “Auto”.
Fast Vol.**	ON	50-950 µL	Activates “Fast vol.” Set the volumes.
	OFF*		Deactivates “Fast vol.”
Blowout	ON		Activates “Repeated blow-out”.
	OFF*		Deactivates “Repeated blow-out”.

\* Default

\*\* The set values correspond to the volume range of the device, e.g. 50 µL-1000 µL.

#### 4.9.2 Parameters in the “Settings”/“Languages” Menu

Parameter	Settings	Explanation
Languages	English* Deutsch Zhōngwén Français Русский	Sets the menu language of the display to the selected language.
* Default		

#### 4.9.3 Parameters in the “Settings”/“Theme” Menu

Parameter	Settings	Explanation
Theme	Default*	Resets the display colour to the defaults.
	Gray, white, green, yellow, orange, red, violet, blue	Sets the colour of the display.
* Default		

#### 4.9.4 Parameters in the “Settings”/“Bluetooth” Menu

Parameter	Settings	Explanation
Bluetooth	ON*	Activates Bluetooth on the device.
	OFF	Deactivates Bluetooth on the device.
Pairing	ON	Activates automatic connection with a known device over Bluetooth.
	OFF*	Deactivates automatic connection with a known device over Bluetooth.
Bluetooth Passkey		The Bluetooth password required to connect is displayed.
* Default		

#### 4.9.5 Parameters in the “Settings”/“Password” Menu

Parameter	Settings	Explanation
Password	ON	Activates password protection.
	OFF*	Deactivates password protection.
Admin	Enter password	Creates the admin password.
User	Enter password	Creates the user password.
* Default		

#### 4.9.6 Parameters in the “Settings”/“Sound” Menu

Parameter	Settings	Explanation
Scrollwheel	ON*	Switches the sound on when the scrollwheel is turned.
	OFF	Switches the sound off when the scrollwheel is turned.
Buttons	ON*	Switches the sound on when a softkey is pressed.
	OFF	Switches the sound off when a softkey is pressed.
Messages	ON*	Switches the sound on when a message is displayed about a low battery level.
	OFF	Switches the sound off when a message is displayed about a low battery level.

\* Default

#### 4.9.7 Parameters in the “Settings”/“Backlight” Menu

Parameter	Settings	Explanation
Backlight	Very Low Low Medium High	Adapts the display lighting based on the parameters selected.

\* Default

#### 4.9.8 Parameters in the “Settings”/“Date & time” Menu

Parameter	Settings	Explanation
Date	DD.MM.YYYY.	Sets the current date.
Time	HH:MM	Sets the current time.

#### 4.9.9 Parameters in the “Settings”/“Tip Ejection” Menu

Parameter	Settings	Explanation
Double Click	ON	Activates ejection of the filter tip by double-clicking on the operating button.
	OFF*	Deactivates ejection of the filter tip by double-clicking on the operating button.
Allow liquid	ON	Activates the ejection of the tip with the safety volume inside.
	OFF*	Deactivates the ejection of the tip with the safety volume inside.

\* Default

#### 4.9.10 Parameters in the “Settings”/“Adjustment” Menu

Parameter	Parameter	Settings	Explanation
Adjustment	Factory		Resets the device adjustment to the default.
	Empty Slot 1	1 point	Set the 1-point adjustment. Set the adjustment point and the adjustment volume. The adjustment volume is freely selectable. The adjustment volume is set on delivery at 10% of the nominal volume.
	Empty Slot 2		
	Empty Slot 3	2 points	Set the 2-point adjustment. Presets the adjustment at 10% and 100% of the nominal volume.
	Empty Slot 4		
Empty Slot 5			
		3 points	Set the 3-point adjustment. Presets the adjustment at 10%, 50% and 100% of the nominal volume.

#### 4.9.11 Parameters in the “Settings”/“User ID” Menu

Parameter	Settings	Explanation
User ID	Enter name:	Creates the user ID. User ID is displayed during device startup.

## 4.9.12 Parameters in the “Settings”/“Reminder” Menu

Parameter	Parameter	Settings	Explanation
Calibration   Maintenance   Quick Check	Last Execution	DD.MM.YY	Displays the date of the last maintenance, calibration or Quick Check (editable).
	Enable	ON	Activates the reminder function for calibration, maintenance, or quick check.
		OFF	Deactivates the reminder function for calibration, maintenance, or quick check..
	Type	Interval	Allows you to enter the period of time until the next calibration, maintenance, or quick check as an interval.
		Date	Allows you to enter a fixed date for the next calibration, maintenance, or quick check.
	Expiration date	DD.MM.YY	Specifies the exact date of the calibration, maintenance, or quick check. Only displayed if [Date] was selected under “Type”.
	Interval value	1 - 4 weeks	Specifies the period of time until the calibration, maintenance, or quick check. Only displayed if [Interval] was selected under “Type”.
		1 - 12 months	
	Snooze	ON	Activates the snooze function for the reminder alarm.
		OFF	Deactivates the snooze function for the reminder alarm.
Lock	Manual	Sets the manual locking function. The device can be manually locked once the date for the next calibration, maintenance, or quick check has arrived.	
	Auto	Sets the automatic locking function. The device is automatically locked once the date for the next calibration, maintenance, or quick check has arrived.	

## 4.9.13 “Settings”/“Reset” Menu

Parameter	Settings	Explanation
Reset	NO	Does <b>not</b> reset the device to the defaults.
	YES	Resets the device to the defaults.

## 5 Installation

### 5.1 Scope of Delivery

Item	Quantity
Picus® 2	1
USB connection cable	1
For models with a volume of > 10 µL: – Safe-Cone Filter – Tweezers	1
For 1-channel models: autoclavable grease	1
ID Stickers	3
Quick Start Guide	1
Quality Control Certificate	1

### 5.2 Unpacking

#### Procedure

- ▶ Unpack the device.
- ▶ Sartorius recommends keeping the original packaging so that the device can be returned properly, e.g. for repairs or maintenance.

### 5.3 Acclimatization

When a cold device is brought into a warmer area: The temperature difference can lead to condensation from humidity in the device (moisture formation). Moisture in the device can lead to malfunctions.

#### Procedure

- ▶ Allow the device to acclimatize to the installation site.

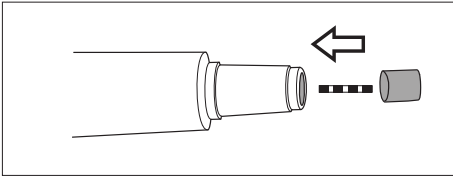
## 5.4 Inserting or Removing a Safe-Cone Filter

### 5.4.1 Inserting a Safe-Cone Filter

Sartorius recommends the use of Safe-Cone filters to avoid contamination. Safe-Cone filters must not be used together with the Safetyspace® filter tips.

#### Procedure

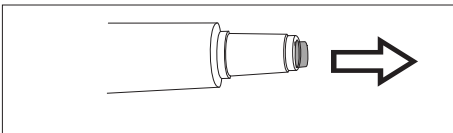
- ▶ Insert a Safe-Cone Filter in the tip cone using the tweezers supplied.



### 5.4.2 Removing the Safe-Cone Filter

#### Procedure

- ▶ Remove the Safe-Cone Filter from the tip cone using the tweezers supplied and dispose of it.





## 6 Commissioning

### 6.1 Clean the Device

It is possible that the device becomes dirty during transport. We recommend cleaning and decontaminating the device before first use (see Chapter "9 Cleaning and Maintenance", page 46).

### 6.2 Charging the Device

The device is delivered with the battery partially charged. We recommend fully charging the device for 1 hour before first use.

#### 6.2.1 Charging with the Charging Carousel or Charging Stand

##### Procedure

- ▶ Make sure the charging stand | charging carousel is connected to the power supply.
- ▶ Insert the device into the charging grooves.
- ▶ Check that the charging contacts of the device are fully inserted in the charging grooves.

#### 6.2.2 Charging with the USB Charger

##### Procedure

- ▶ Slide open the cover of the USB port (1).
- ▶ Connect the connection cable to the micro USB port of the device.
- ▶ Connect the other end of the connection cable to a mains plug via the USB- port
- ▶ Connect the mains plug to the socket (supply voltage) at the installation site.



## 6.3 Switching the Device On or Off

### Procedure

- ▶ To switch the device on:
  - ▶ Press the On | Off button.
  - ▶ If password protection is activated: Log in with the password.
  - ▶ When the device is switched on for the first time: Enter the date and the time in the query.
- ▷ The device actuates the piston.
- ▷ The device is ready for use.
  
- ▶ To switch the device off:
  - ▶ If password protection is activated: Log out of the profile being used. To do this, briefly press the On | Off button.
  - ▶ Press and hold the On | Off button for a few seconds.

### 6.3.1 Returning to Active Mode

During use and charging, the device is in active mode. All processor functions are activated and the backlighting on the display is switched on. When the device is not used (for a long time) it behaves as follows:

- If the device is not used for more than 30 seconds: The device is in energy saving mode and the backlight is dimmed. Pressing any key or turning the scrollwheel let return to the active mode.
- If the device is not used for more than 10 minutes: The backlight switches off. Pressing the On | Off button lets return to active mode.
- If the device is not used for more than 4 hours: The device switches itself off.

## 6.4 Rinsing Tips

Sartorius recommends to rinse the tips before operating.

### Requirements

Tips are attached to the cone.

### Procedure

- ▶ Open the main menu.
- ▶ Set the operating modus "Pipetting".
- ▶ Aspirate water. To do so, press the operating button.
- ▶ Dispense water. To do so, press the operating button again.
- ▶ Repeat aspirating and dispensing 3 - 5 times.
- ▷ The Tips are rinsed.

# 7 System Settings

## 7.1 Setting Defaults

Defaults, which are coordinated with the ambient conditions and the requirements during operation, can be set for the device and the applications.

To configure the device, the following settings are recommended:

- Set the menu language
- Set the adjustment

### Procedure

- ▶ Open the main menu.
- ▶ Open the “Settings” sub-menu.
- ▶ To make settings: Open the desired menu item.
- ▶ Select the desired parameters and confirm (for parameters, see Chapter “4.9 List of Parameters”, page 26).
- ▶ Exit the menu.

## 7.2 Using the Device with the Sartorius Pipetting MobileApp

### 7.2.1 Downloading the Sartorius Pipetting MobileApp

#### Procedure

- ▶ Download the Sartorius Pipetting MobileApp. To do this, select one of the following options:
  - ▶ Scan the QR code on the Quick Start Guide.
  - ▶ Open the App Store on the device and find the Sartorius Pipetting MobileApp.

### 7.2.2 Adding the Device in the app

#### Requirements

Bluetooth® is activated on the device (see Chapter “4.9.4 Parameters in the “Settings”/“Bluetooth” Menu”, page 27).

- ▶ When the app has been successfully downloaded: Open the Sartorius mobile app.
- ▶ Create Sartorius ID on <https://my.sartorius.com/>.
- ▶ Enter the Sartorius ID and the password.
- ▷ The main menu is displayed.
- ▶ Open the “PIPETTE MANAGEMENT” menu.
- ▶ Add the desired device. To do this, proceed as follows:
  - ▶ Press the [+] button.
  - ▶ Follow the wizard’s instructions.
- ▷ The device is listed in the “List of all pipettes” menu.

## 7.2.3 Creating the Pipette Set

### Requirement

Devices are added in the Sartorius Pipetting MobileApp (see Chapter “7.2.2 Adding the Device in the app”, page 35).

### Procedure

- ▶ Open the “MY PIPETTE SETS” menu.
- ▶ Press the [+] button.
- ▷ The “New pipette set” display appears.
- ▶ Follow the wizard’s instructions.
- ▷ The pipette set is listed in the “MY PIPETTE SETS” menu.

## 7.2.4 Starting the Workflow

### Requirement

- Devices are added in the app (see Chapter “7.2.2 Adding the Device in the app”, page 35).
- Bluetooth® is activated on the device (see Chapter “4.9.4 Parameters in the “Settings”/“Bluetooth” Menu”, page 27).

### Procedure

- ▶ Open the “WORKFLOWS” menu.
- ▶ Select the desired workflow.
- ▶ Follow the wizard’s instructions.

## 7.3 Performing a Software Update

A software update can be performed via the Sartorius Pipetting mobile app. In the Sartorius Pipetting mobile app, you can see which devices have not yet been updated to the current software version.

The current software version of the device is specified under the parameter “Settings”/“Information”.

A software update can extend or change the functionality of the device. Sartorius recommends performing the software update regularly.

### Requirement

- The desired device is switched on.
- If password protection is activated: Be logged in with the admin password.
- Bluetooth is activated on the device (see Chapter “4.9.4 Parameters in the “Settings”/“Bluetooth” Menu”, page 27).
- The device is added to the Sartorius Pipetting Mobile app (see Chapter “7.2.2 Adding the Device in the app”, page 35).

### Procedure

- ▶ Open the Sartorius Pipetting mobile app and log in with your Sartorius ID.
- ▶ Open the menu "PIPETTE MANAGEMENT".
- ▷ The App searches for devices nearby.
- ▷ A list of found devices opens.
- ▶ Select the desired device and tap on the [UPDATE] button.
- ▷ The device display displays the update progress.
- ▷ When the update is completed, the [UPDATE] button is no longer displayed.

## 7.4 Managing User Data

### 7.4.1 Creating a User ID

#### Procedure

- ▶ Open the main menu.
- ▶ Open the "Settings" sub-menu.
- ▶ Open the "User ID" menu option.
- ▶ Create a name and confirm.

### 7.4.2 Activating Password Protection

The device has two levels of password protection. The top level is for the administrator with all access rights. The bottom level is for the user with limited access rights. If password protection is active: Only the administrator has access to the following functions:

- Managing Bluetooth® settings
- Managing passwords
- Setting the date and time
- Setting reminders
- Creating and saving protocols
- Saving programs (hotkey)
- Managing adjustment settings
- Resetting to factory default settings

When the password is activated, users have limited access to the device functions and can call up the following functions, for example: operating modes, backlight or user ID.

#### Procedure

- ▶ Open the main menu.
- ▶ Open the "Settings" sub-menu.
- ▶ Open the "Password" menu option.
- ▶ Select the "Login" parameter and set to "ON".
- ▶ Select the parameters "User" and "Administrator" one after the other and assign a password.
- ▶ Confirm the entry with the [OK] soft key.
- ▶ Save the changes made with the [SAVE] soft key.
- ▷ Password protection is enabled.

### 7.4.3 Logging Into or Out of the Device

When password protection is activated, operators must log into and out of the device. Login is performed directly after switching on. Logging out is **not** possible during an active operating mode.

#### Procedure

- ▶ To log in the operator: Switch on the device.
- ▷ A password request is displayed.
- ▶ Enter the user or administrator password and confirm.
- ▶ To log out the operator: Briefly press the On | Off button.
- ▷ The user profile is logged out.

### 7.4.4 Deactivating Password Protection

#### Requirements

The administrator profile is logged in.

#### Procedure

- ▶ Open the main menu.
- ▶ Open the "Settings" sub-menu.
- ▶ Open the "Password" menu option.
- ▷ A password request appears.
- ▶ Enter the administrator password.
- ▶ Select the parameter "Password" and set to "OFF".
- ▶ Save the changes made with the [SAVE] soft key.
- ▷ Password protection is deactivated.

## 7.5 Adjusting the Device

The piston stroke is designed for normal conditions as standard:

- Aqueous liquids
- Normal pressure
- Device, tip and liquid at room temperature

If the conditions change: The accuracy of the dispensed volume may change. The adjustment function can ensure that the device remains accurate. Adjustment of the dispensed volume is required in the following cases:

- The pipetting liquid's properties are significantly different from those of water, e.g. viscous or volatile liquids.
- The temperature of the device, pipetting tip and liquid deviates significantly.
- The ambient air pressure differs from normal pressure.
- The routine test shows a measurement deviation from the test specifications of the pipette (routine test see Chapter "9.6 Performing the Routine Test", page 56).

The adjustment function enables the device to be set to one or more adjustment points. The more adjustment points are selected, the greater the accuracy over the entire volume range of the device. Sartorius recommends the following settings:

- 1-point adjustment: Pipetting a constant volume within a range
- 2-point or 3-point adjustment: Pipetting volumes within the entire range

When adjusting the accuracy, the actual volumes obtained should be measured in pipetting mode. After adjustment has been carried out, it will be applied to all modes and a symbol for adjustment will be displayed.

### Procedure

- ▶ If password protection is activated: Log in with the admin password.
- ▶ Open the main menu.
- ▶ Open the "Settings" sub-menu.
- ▶ Open the "Adjustment" menu option.
- ▶ Select the parameter under which the adjustment settings should be saved, e.g. Empty Slot 1.
- ▶ Activate edit mode.
- ▶ Set a name.
- ▶ Select the number of adjustment points.
- ▶ Confirm the selection.

### Set the target volume for adjustment points (1-point adjustment)

#### Procedure

- ▶ For 1-point adjustment: Set the target volume and press [EDIT].



### Set the target volume for adjustment points (2-point or 3-point adjustment)

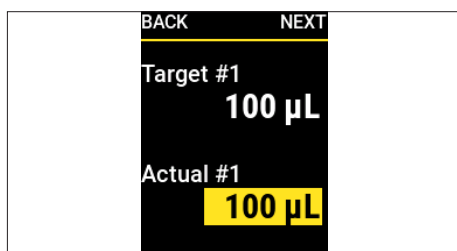
#### Procedure

- ▶ Confirm the automatically set target volume.

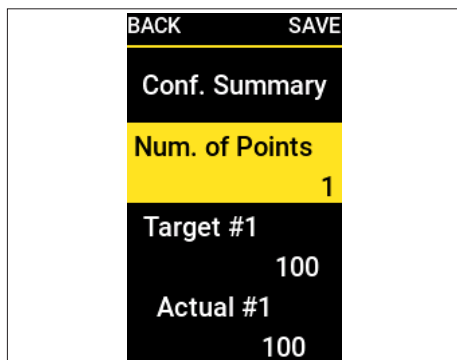
### Setting the measured volume

#### Procedure

- ▶ Enter the volume actually measured and confirm.



- ▶ Configuration summary is displayed.
- ▶ Save the set adjustment data.

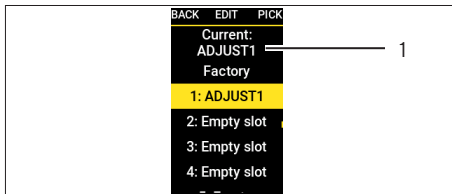




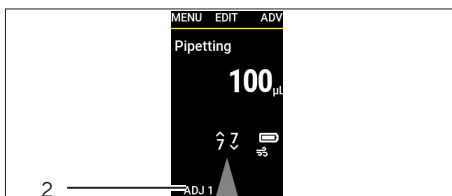
## 7.5.1 Activating the Adjustment Setting

### Procedure

- ▶ If password protection is activated: Log in with the admin password.
- ▶ Open the main menu.
- ▶ Open the “Settings” sub-menu.
- ▶ Open the “Adjustment” menu option.
- ▶ Pick the desired parameter.
- ▷ The top of the menu (1) indicates which parameter is activated.



- ▶ Return to operating mode.
- ▷ The adjustment is applied to all operating modes.
- ▷ The selected adjustment setting is displayed (2).



## 7.5.2 Documenting Adjustment for Liquids Other Than Water

### Procedure

- ▶ If the device was adjusted for use with liquids other than water: Attach a label to the outside of the device indicating the name of the liquid and the adjusted volume range. To do this, the supplied ID stickers can be used.

## 7.6 Setting Up Reminders

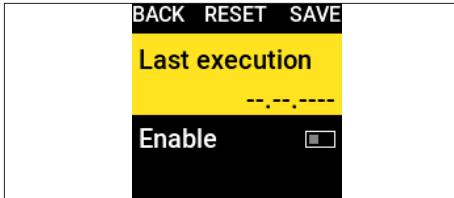
The device enables the configuration of reminders for the following activities:

- Maintenance
- Calibration
- Quick checks

Reminders can be set as a date or interval (weeks |months). To set both a date and an interval: The criterion that is reached first triggers the reminder function and reminds the user about the due measures via a message and a signal sound. It is possible to snooze the reminder for a set number of days. For deadline-based reminders, the user is reminded 14 days before the deadline.

### Procedure

- ▶ If password protection is activated: Log in with the admin password.
- ▶ Open the main menu.
- ▶ Open the “Settings” sub-menu.
- ▶ Open the “Reminder” menu option.
- ▶ Select and confirm the reminder type, e.g. calibration.
- ▶ Set the reminder functions one after the other:
  - ▶ Set the date of the last calibration.
  - ▶ If necessary: Define the date of the next calibration and apply.
  - ▶ If necessary: Define the interval until the next calibration and apply.
  - ▶ If necessary: Activate the snooze function and set the number of days.
- ▶ When all desired reminder functions have been set: Press the [SAVE] softkey.
- ▶ All settings are saved and the reminder function is activated.
- ▶ The display returns to the “Settings” sub-menu.



## 7.6.1 Deactivating Reminders

### Procedure

- ▶ If password protection is activated: Log in with the admin password.
- ▶ Open the main menu.
- ▶ Open the “Settings” sub-menu.
- ▶ Open the “Reminder” menu option.
- ▶ Select and confirm the reminder type.
- ▶ To deactivate a reminder function: Set to [OFF] and skip to the next parameter.
- ▶ When the desired reminder functions have been deactivated: Press the [SAVE] softkey.
- ▶ All settings are saved and the reminder function is deactivated.
- ▶ The display returns to the “Settings” sub-menu.

# 8 Operation

## 8.1 Accessing Digital Compliance Labels

Digital compliance labels can be displayed in the device. These show which country-specific regulations and standards the device complies with. More detailed information on the conformity of the device can be found in the chapter “Conformity of the Device” (see Chapter 17, page 76).

### Procedure

- ▶ Open the main menu.
- ▶ Open the menu entry “Compliance”.
- ▷ The information on conformity including the digital labels is displayed.
- ▶ If necessary: Scroll down to see all labels.

## 8.2 Using Operating Modes (Examples)

### 8.2.1 Multi Dispensing

#### Requirements

Tips are attached to the cone.

#### Procedure

- ▶ Open the main menu.
- ▶ Select the “Multi Disp.” pipetting mode and confirm.
- ▷ The display shows the last pipetting settings for the “Multi Disp.” pipetting mode.
- ▶ To change pipetting settings: Activate edit mode.
- ▶ Edit the desired settings and confirm.
- ▶ To activate advanced functions or adapt the safety volume:
  - ▶ Press the [ADV] softkey.
  - ▶ Select and activate the desired advanced functions (see Chapter “4.7 Menu Structure in the “Advanced Functions” Menu”, page 25).
- ▶ To aspirate liquid: Press the operating button.
- ▷ The selected volume and safety volume are aspirated.
- ▶ To eject the primary excess: Press the operating button.
- ▶ Press the operating button until all sub-volumes have been dispensed.
- ▷ The message: “Double-click to empty?” is displayed.
- ▶ To empty the tip: Press the operating button twice.
- ▶ To eject the tip: Press the electronic tip ejection.

## 8.2.2 Manual Pipetting

### Requirements

Tips are attached to the cone.

### Procedure

- ▶ Open the main menu.
- ▶ Select the “Manual” pipetting mode.
- ▶ To change pipetting settings: Activate edit mode.
- ▶ Edit the desired settings and confirm.
- ▶ To start the pipetting process: Press the operating button.
- ▶ Aspirate the liquid. To do this, use one of the following options:
  - ▶ Press and hold the operating button. The aspiration speed is constant.
  - ▶ Turn the scrollwheel counterclockwise. The further the scrollwheel is turned, the higher the aspiration speed.
- ▶ To interrupt the process: Release the operating button or the scrollwheel.
- ▶ To switch to dispensing: Turn the scrollwheel briefly clockwise.
- ▷ The display arrow points down.

### Dispensing liquids

#### Procedure

- ▶ Dispense the liquid. To do this, use one of the following options:
  - ▶ Press and hold the operating button. The dispensing speed is constant.
  - ▶ Turn the scrollwheel clockwise. The further the scrollwheel is turned, the higher the dispensing speed.
- ▶ When the entire volume has been dispensed: Press the operating button.
- ▷ The message “Double-click to empty?” is displayed.
- ▶ To empty the tip: Press the operating button.
- ▶ To eject the tip: Press the electronic tip ejection.

## 8.3 Locking and Unlocking the Device

### 8.3.1 Activating the Device Lock

The device can be locked to prevent further use. This function can be used when an error has occurred in the device, for example, or the device has become contaminated.

#### Procedure

- ▶ Open the main menu.
- ▶ Open the "Settings" sub-menu.
- ▶ Select the "Lock" menu item and confirm.
- ▷ The display shows "Lock pipette"?
- ▶ Confirm the lock with the [YES] softkey.
- ▷ The device is locked.

### 8.3.2 Opening the Device Lock

#### Procedure

- ▶ Press the [UNLOCK] softkey.
- ▶ If password protection is activated: Enter the admin password.
- ▷ The device is unlocked and the device is ready for use.

## 9 Cleaning and Maintenance

### 9.1 Cleaning the Surfaces of the Device

We recommend to clean the surface of the device daily.

- Materials:
- Gloves
  - Cleaning agents
  - Soft, lint-free cleaning cloth

#### Requirements

- The process is finished.
- The filter tip has been ejected.

#### Procedure

- ▶ Switch off the device.
- ▶ Only use suitable cleaning agents and cleaning procedures and note the product information for the cleaning agent used (see Chapter “14.10.1 Cleaning Agents, Decontamination Agents and Cleaning Methods”, page 68).
- ▶ Dampen the cleaning cloth with the cleaning agent.
- ▶ Clean the device housing with the damp cleaning cloth.
- ▶ Wipe the device housing dry.

### 9.2 Maintenance Schedule

Interval	Component	Activity	Chapter, page
Regularly, depending on the operating conditions	Safe-Cone Filter	Replace the Safe-Cone Filters.	5.4.2, 32
	1-channel models: Lower assembly of the device	Clean the lower assembly and grease components.	9.3, 47
		Sterilise the lower assembly	9.4, 52
Multi-channel models: Lower assembly of the device	Sterilise the lower assembly	9.4, 52	
	For cleaning and greasing the lower assembly contact Sartorius Service.		
Regularly, e.g., every 3 months and after each in-house maintenance	Device	Perform a routine test to check the pipette performance.	9.6, 56
6 to 24 months, depending on the operating conditions	Device	Contact Sartorius Service for calibrating the device.	

## 9.3 Cleaning and Greasing the Lower Assembly of the Device

### 9.3.1 Preparing the Device

The working steps described for cleaning and greasing the lower assembly only apply to 1-channel models. For cleaning and greasing multi-channel pipettes contact Sartorius Service.

#### Procedure

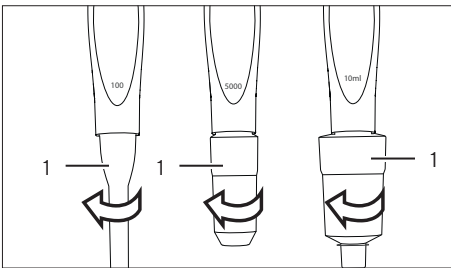
- ▶ **⚠ CAUTION** Risk of injury due to unprotected, moving parts! Exposed moving parts can cause injuries. Switch off the device before cleaning and maintenance work.
- ▶ Remove the Safe-Cone Filters (see Chapter 5.4.2, page 32).

### 9.3.2 Dismantling the Lower Assembly (1-channel Models)

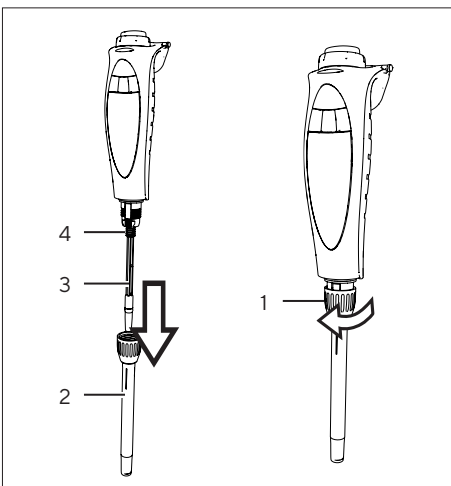
#### Models with maximum volume of up to 1,000 $\mu\text{L}$

#### Procedure

- ▶ Unscrew the tip ejector (1) in a counterclockwise direction.



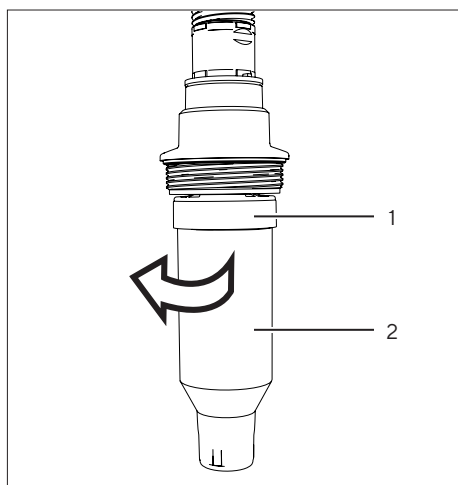
- ▶ Turn the tip cone holder (1) counterclockwise and slowly pull off together with the tip cone (2).
- ▶ Slowly pull the spring (4) off the piston (3).



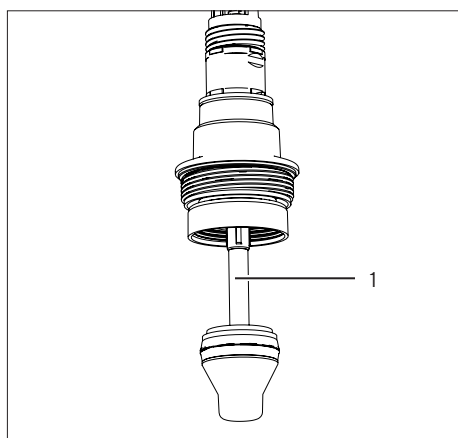
Models with maximum volume of 5,000  $\mu$ L

Procedure

- ▶ Hold the tip cone holder (1) with one hand and use the other hand to turn the tip cone cylinder (2) counter-clockwise and pull it off.



- ▶ The piston (1) is exposed.







### 9.3.4 Greasing Components (1-channel models)

---

Materials:     – Gloves  
                   – Grease (included in scope of delivery)

---

Tool:            Brush

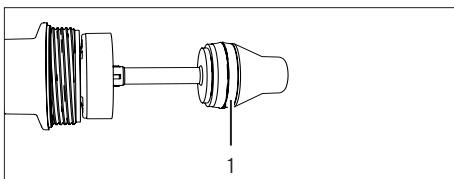
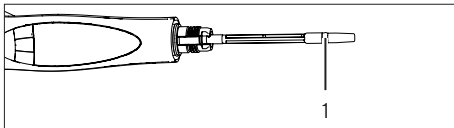
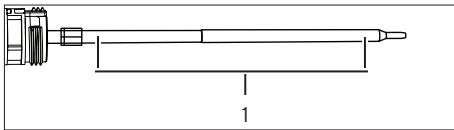
---

#### Requirements

The lower assembly is dismantled.

#### Procedure

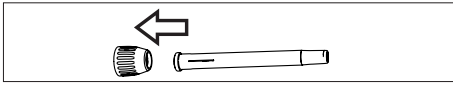
- ▶ Wet the brush with the grease supplied. To do this, proceed as follows:
  - ▶ Models with maximum volume of 10  $\mu\text{L}$  | 120  $\mu\text{L}$ : Apply a thin layer of grease to the piston (1).
  - ▶ Models with maximum volume of 300  $\mu\text{L}$  | 1,000  $\mu\text{L}$ : Apply a thin layer of grease around the seal (1).
  - ▶ Models with maximum volume of 5,000  $\mu\text{L}$  | 10,000  $\mu\text{L}$ : Apply a thin layer of grease inside the tip cone and around the seal (1).



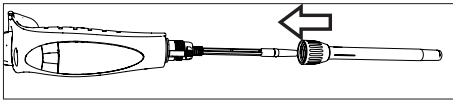
### 9.3.5 Assembling the Lower Assembly (1-channel Models)

#### Models with maximum volume of up to 1,000 $\mu\text{L}$

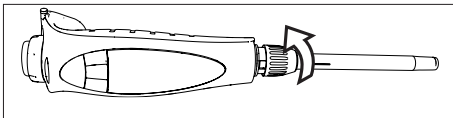
##### Procedure



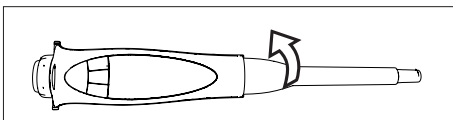
- ▶ Slide the spring over the piston.
- ▶ Place the tip cone in the tip cone holder.



- ▶ Slide the tip cone holder together with the tip cone over the piston.



- ▶ To secure the tip cone holder and tip cone: Turn the tip cone holder clockwise to tighten.

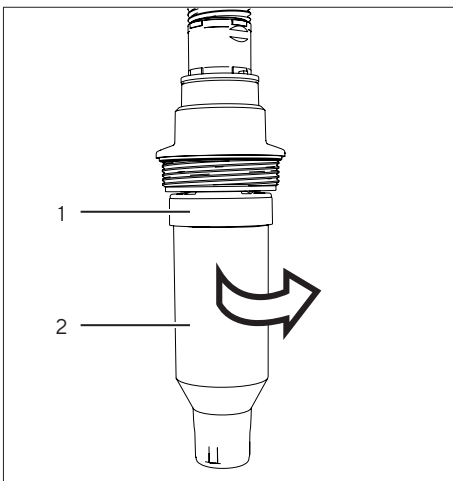


- ▶ Slide the tip ejector over the tip cone and secure by turning it clockwise.

- ▶ Insert a new Safe-Cone Filter (see Chapter 5.4, page 32).
- ▶ To ensure that the grease is evenly distributed:
  - ▶ Switch on the device.
  - ▶ Press the operating button several times.
- ▶ Check the operation of the device.

#### Models with maximum volume of 5,000 $\mu\text{L}$

##### Procedure

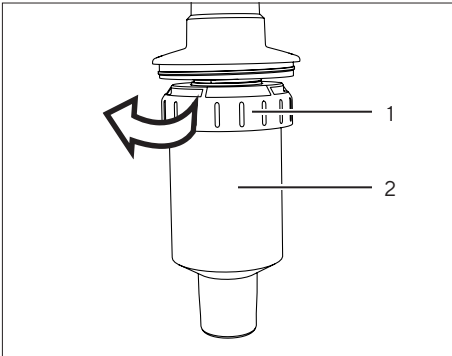


- ▶ To secure the tip cone: Hold the tip cone holder (1) with one hand and use the other hand to turn the tip cone cylinder (2) clockwise.
- ▶ Check that the tip cone is **not** too tight.

## Models with maximum volume of 10,000 µL

### Procedure

- ▶ Slide the tip cone cylinder (2) slowly over the piston.
- ▶ Turn the locking ring (1) counterclockwise.
- ▶ Check that the tip cone is **not** too tight.



## 9.4 Sterilising the Device

### 9.4.1 Selecting the Sterilization Method

#### Procedure

- ▶ Select the preferred sterilization method for the device (see Chapter “14.10.2 Approved Sterilization Methods for Lower Assembly”, page 68).

### 9.4.2 Sterilizing the Device with UV Radiation

The device is made of UV-resistant materials and tolerates temporary exposure to UV radiation. A prolonged or frequent exposure to UV radiation may cause yellowing and brittling of the device.

#### Procedure

- ▶ Sterilize the lower assembly with UV radiation.

### 9.4.3 Sterilizing the Device with a Decontamination Agent

---

Materials:      – Gloves  
                      – Suitable decontamination agent (see Chapter 14.10.1, page 68)

---

Tool:             Soft, lint-free cleaning cloth

---

#### Procedure

- ▶ Dampen the cleaning cloth with the decontamination agent.
- ▶ During decontamination work, ensure that **no** liquids get inside the device.
- ▶ Wipe the outside of the device with the damp cleaning cloth and leave it to dry.

## 9.4.4 Sterilising the Lower Assembly of the Device

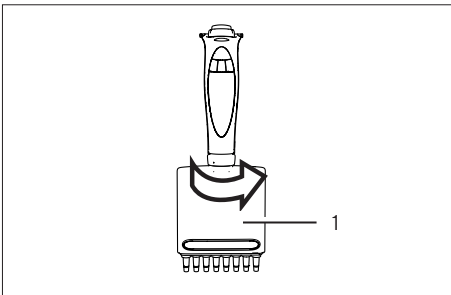
Materials: Autoclaving bag

### Requirements

The device is autoclavable (see Chapter “14.10.1 Cleaning Agents, Decontamination Agents and Cleaning Methods”, page 68).

### Procedure

- ▶ Remove the Safe-Cone Filters (see Chapter 5.4.2, page 32).
- ▶ If a 1-channel model is used: Dismantle the lower assembly (see Chapter 9.3.2, page 47).
- ▶ If a multi-channel model is used: Unscrew the lower assembly (1) clockwise and remove it.

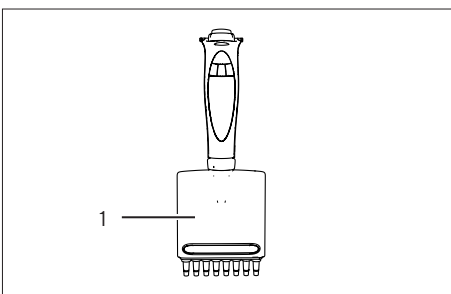


- ▶ Place the components of the lower assembly in the autoclaving bag:
  - 1-channel model: Tip ejector, tip cone, tip cone holder
  - Multit-channel model: Lower assembly
  - Autoclave the components in the autoclaving bag. Comply with the autoclaving conditions (see Chapter “14.10.2 Approved Sterilization Methods for Lower Assembly”, page 68).
- ▶ Leave the components to cool and dry.

### Finishing the Sterilization by Autoclaving

#### Procedure

- ▶ Assemble all components on the lower assembly. To do that proceed as follows:
  - ▶ If a 1-channel model is used: Assemble the lower assembly (see Chapter 9.3.5, page 51).
  - ▶ If a multi-channel model is used: Screw the lower assembly (1) onto the control head by turning it counterclockwise.



## 9.5 Defining the Test Routine for the Routine Test

We recommend testing the performance of the pipette regularly in a routine test (for intervals, see Chapter “9.2 Maintenance Schedule”, page 46).

We recommend defining a test routine for the routine test that takes into account the following criteria:

Test Routine	Criteria	Explanation
General test routine	Accuracy requirements of the relevant application	
	Frequency of use	
	Number of operators for the device	
	Nature of the liquid dispensed	
	Maximum permitted error limits for systematic and random errors, as per ISO 8655-2	As per ISO 8655-2
	Acceptable error range taking into account: <ul style="list-style-type: none"> <li>– Application</li> <li>– Application area</li> <li>– Conditions for precision of the application</li> </ul>	Sartorius specifications were achieved in strictly controlled conditions, as per ISO 8655-6 (see Chapter “14.13 Performance Specification”, page 70).
	Multistage routine test with three volumes recommended: <ul style="list-style-type: none"> <li>– 100% of the nominal volume</li> <li>– 50% of the nominal volume</li> <li>– 10% of the nominal volume</li> </ul>	
Risk of the application, e.g. in relation to accuracy		
Additional requirements for multi-channel pipettes	All channels of the multi-channel pipette must be tested individually. The following options are available for doing this:	
	Use a multi-channel balance that measures the dispensed liquid from all channels in parallel.	The test liquid is aspirated, dispensed and measured in all channels at the same time.
	Use a single-channel balance that measures the dispensed test liquid from one channel.	The test liquid for each channel is aspirated, dispensed and measured individually in succession. The test liquid from the other channels is discarded as residual water.

Test Routine	Criteria	Explanation
Additional requirements for testing in the multi dispensing pipetting mode	Required test volume	Nominal volume, with 10 measurements with 10% of the nominal volume
	Excess liquid volume	Must be discarded after aspirating
	Pause after dispensing of the first liquid volume	The test liquid does <b>not</b> move during this process.
	Changing the tip depending on the amount of test liquid	<ul style="list-style-type: none"> <li>– If there is enough test liquid for 10 measurements: Do <b>not</b> change the tip.</li> <li>– If there is <b>not</b> enough test liquid for 10 measurements: <ul style="list-style-type: none"> <li>– change the tip between the measurements.</li> <li>– After changing, once again dispense excess liquid volume, which is discarded.</li> </ul> </li> </ul>
	Residual liquid	Must be discarded.

#### Procedure

- ▶ Prepare a test routine that takes into account the required criteria for the routine test on the device.

## 9.6 Performing the Routine Test

A routine test with an analytical balance (single-channel balance) is described below. Further information on the possible test routines is described in the ISO 8655 series of standards, e.g., for the testing of multi dispensing in pipetting mode.

- Materials:
- Test liquid (test water): distilled, deionized water (ISO 3696, quality 3)
  - Test water container
  - Measuring container, for positioning on the analytical balance
  - Pipette tips
- Tool: Analytical balance, in compliance with the standard ISO 8655-6

### Requirements

The test water container is filled with test water.

### Procedure

- ▶ Check whether the following conditions are met:
  - The required ambient conditions have been complied with (see Chapter “14.2 Ambient Conditions”, page 64).
  - The device, tips and test water have been stored in the room for at least two hours so that they have adjusted to the ambient conditions.
  - The “Pipetting” mode is set in the device.
  - The aspirating speed and dispensing speed are set to 7.
- ▶ If necessary: Ensure that the conditions are complied with or adjust the settings on the device.

### 9.6.1 Aspirating the Test Water

The conditions for aspirating the test water must be complied with (see Chapter “14.12.1 Conditions for Aspirating the Test Liquid”, page 69).

### Procedure

- ▶ Adjust the preferred test volume ( $V_S$ ).
- ▶ Apply the tip to the tip cone.
- ▶ To achieve a humidity balance in device: Aspirate and dispense the test water 5 times with the tip.
- ▶ Aspirate the test water. To do this, proceed as follows:
  - ▶ Hold the device vertically, press the control button and immerse the tip below the surface of the water to the specified immersion depth.
  - ▶ Comply with the specified wait time.
- ▶ Remove the device from the test water container vertically.



## 9.6.2 Dispensing Test Water and Performing the Measurement

We recommend performing 10 measurements. The tip must be replaced after 5 measurements.

### Requirements

The analytical balance is ready for measuring.

### Procedure

- ▶ Dispense the test water. To do this, proceed as follows:
  - ▶ Hold the tip above the liquid level at an angle of 30° - 40° on the inner wall of the measuring container.
  - ▶ Press and hold the control button and release the test water into the measuring container.
  - ▶ In order to remove drops on the tip: Drag the tip 8 - 10 mm along the inner wall of the measuring container.
- ▶ Release the control button.
- ▶ Remove the device from the measuring container vertically.
- ▶ Read off the weight in mg ( $m_i$ ).
- ▶ Repeat the test cycle until 10 measurements have been recorded. Replace the tip after five test cycles.

## 9.6.3 Evaluating Measured Values

### Requirements

Ten measurements have been recorded.

### Procedure

- ▶ Convert the recorded masses ( $m_i$ ) into volume ( $V_i$ ) by multiplying with a correction factor  $Z$ . The following applies here:  $V_i = m_i \cdot Z$  (for correction factor  $Z$ , see Chapter 14.12.2, page 69).
- ▶ Calculate the average volume: ( $V$ ):  $V = \Sigma(V_i)/10$ .
- ▶ To evaluate conformity: Calculate the systematic error " $e_s$ " of the measurement. Use the following formula to do this:
  - In  $\mu\text{L}$ :  $e_s = V - V_s$
  - Or in %:  $e_s = 100 (V - V_s)/V_s$
  - $V_s$  = selected test volume
- ▶ For the conformity assessment, calculate the random measurement deviation as standard deviation or as variation coefficient:
  - As standard deviation ( $n$  = number of measurements (10))
  - As variation coefficient  $C_v = 100\% \text{ sr} / V$

$$s = \sqrt{\frac{\Sigma(V_i - \bar{V})^2}{n - 1}}$$

- ▶ Compare the systematic error and the random error with the values of the performance specification from your own laboratory.
- ▷ If the measurement results are within the performance specification: The device is ready for use.
- ▷ If the measurement results are **not** within the performance specification:
  - ▶ Check the device for both systematic errors and random errors.
  - ▶ Ensure that a suitable tip is used.
  - ▶ Ensure that a good seal has been formed between tip and tip cone.
  - ▷ If necessary: Adjust the device (see Chapter 7.5, page 39).

# 10 Trouble

## 10.1 Troubleshooting

Fault	Cause	Remedy	Chapter, page
Droplets remain in the tip.	The tip is not compatible.	Use original Sartorius tips. Check that the tip is securely seated and, if necessary, tighten.	
	The device is contaminated.	Clean the device.	9, 46
	The device is defective.	Replace the defective parts. Send the device in for servicing.	
Inaccurate function	The device is contaminated.	Clean the device.	9, 46
	The device is defective.	Replace the defective parts. Send the device in for servicing.	
The device is switched off.	Pipette in Off mode	Switch on the device.	6.3.1, 34
	The battery is empty.	Charge the device.	6.2, 33
The piston jams	Internal components are loose	Open the lower assembly and ensure the parts are securely fitted.	9.3.2, 47
	Pipette contaminated	Clean the pipette.	9, 46
	Pipette defective	Replace defective parts. Send the device in for servicing.	
The aspirated volume is too small	The Safe-Cone Filter is contaminated	Replace the Safe-Cone Filter.	5.4, 32
	The device is contaminated.	Clean the device.	9, 46
	The device is defective	Replace the defective parts. Send the device in for servicing.	
The device does not respond.		Reset the device to the defaults.	10.2, 60
		Send the device in for servicing.	

## 10.2 Resetting to Defaults

When resetting to defaults, the stored settings and the other information contained in the memory are not changed.

### Requirements

- The device is not connected to the charger.
- The device is not connected to the USB charger.

### Procedure

- ▶ Press and hold the On | Off button and the right softkey button simultaneously for a few seconds.
- ▷ The device switches itself off.
- ▶ To switch on: Press the On | Off button.
- ▷ The device is reset to the defaults.

# 11 Decommissioning

## 11.1 Decommissioning the Device

### Requirements

The filter tip(s) has | have been ejected.

### Procedure

- ▶ If password protection is activated: Log out of the user profile.
- ▶ Switch off the device.
- ▶ Remove the Safe-Cone Filter.
- ▶ Clean the device.

# 12 Storage and Shipping

## 12.1 Storage

### Procedure

- ▶ Switch off the device.
- ▶ Check that the ambient conditions are met (see Chapter “14.2 Ambient Conditions”, page 64).
- ▶ Store the device in vertical position, e.g., in a pipette stand or charging stand.
- ▶ If a charging stand or charging carousel is used:
  - ▶ Place the device in the charging stand.
  - ▶ If the device is not used for several months: Disconnect the charging stand from the power supply.

## 12.2 Returning the Device and Parts

Defective devices or parts can be returned to Sartorius. Returned devices must be clean, decontaminated, and packaged properly.

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

Devices contaminated with hazardous substances, e.g. harmful biological or chemical substances, will **not** be accepted for repair and disposal.

### Procedure

- ▶ Decommission the device.
- ▶ If necessary: Decontaminate the device.
- ▶ Contact Sartorius Service for information on how to return devices or parts (see [www.sartorius.com](http://www.sartorius.com)).
- ▶ Pack the device and its parts properly for return.

# 13 Disposal

## 13.1 Disposing of the Device and Parts

The device and its accessories must be disposed of properly by disposal facilities.

A lithium battery is installed inside the device. Batteries must be disposed of properly by disposal facilities.

### Procedure

- ▶ Dispose of the device in accordance with the country-specific legal provisions. Inform the disposal facility that there is a lithium battery inside the device.
- ▶ Dispose of the packaging in accordance with the country-specific legal provisions.
- ▶ Dispose of the consumables in accordance with the country-specific legal provisions.

# 14 Technical Data

## 14.1 Dimensions and Weights

### 14.1.1 1-channel models

		LH-747021	LH-747041	LH-747061	LH-747081	LH-747101	LH-747111
	Unit	Value	Value	Value	Value	Value	Value
Length	mm	208	214	211	214	185	185
Weight	g	102	104	103	105	117	127

### 14.1.2 Multi-channel models with 8

		LH-747321	LH-747341	LH-747361	LH-747391
	Unit	Value	Value	Value	Value
Length	mm	217	218	214	214
Weight	g	167	169	164	176

### 14.1.3 Multi-channel models with 12

		LH-747421	LH-747441	LH-747461	LH-747491
	Unit	Value	Value	Value	Value
Length	mm	217	218	214	214
Weight	g	195	197	190	215

## 14.2 Ambient Conditions

	Unit	Value
Laboratory, for indoor use only		
Pollution level according to IEC 61010-1		II
Maximal height above sea level	m	2000
Suitable for protection class		
Protection class of the device, as per DIN EN 60529-1		IP 40
Temperature		
In operation	°C	+15 - +30
During routine test	°C	+15 - +30
In storage   during transport	°C	-20 - +40
Relative humidity, in operation		
At temperatures up to 31°C, maximum	%	80
<b>No</b> potentially explosive areas		
Other properties		
Store dry		

## 14.3 Electrical Data

### 14.3.1 Power Supply

	Unit	Value
Alternating voltage	V <sub>AC</sub>	100 - 240 (± 10 %)
Frequency	Hz	50 - 60
Current consumption, maximum	A	0.125
Power supply only permitted through mains connection cable or storage devices provided by Sartorius		
Protection class, as per IEC60950-1, (reinforced insulation between mains and secondary circuit)		II
Surge category according to IEC 60664-1		II
Connection cable		
Connection cable with micro-USB port		
Length	m	1,8



### 14.3.2 Safety of Electrical Equipment

---

Safety provisions in accordance with EN 61010-1 / IEC 61010-1 Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements

---

Safety of electrical equipment, in accordance with EN 61326-1 / IEC 61326-1 Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1: General requirements.

---

### 14.3.3 Safety and Electromagnetic Compatibility

---

Safety of electrical equipment, in accordance with EN 61326-1 / IEC 61326-1 Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1: General requirements.

---

Interference immunity: Suitable for use in industrial areas

---

Emitted interference: Class B; Suitable for use in residential areas and areas that are directly connected to a low-voltage network that also supplies residential buildings.

---

## 14.4 Interfaces

	Unit	Value
Micro USB port		
Communication: USB host (Connection cable)		
Bluetooth® 5.3 LE, BMD-350		
Transmission power, maximum	dBm	4
Range	m	10

## 14.5 Operating Modes with Advanced Functions

	Tracker	Mixing	Counter	Safety Volume	Auto-Dispensing	Fast Dispensing	Repeated Blow-out	Auto pre-out
Pipetting	▪	▪	▪				▪	
Reverse Pipetting	▪		▪	▪				
Manual Pipetting							▪	
Multi Dispensing	▪			▪	▪			▪
Diluting		▪					▪	
Sequential Dispensing	▪			▪	▪			▪
Multi Aspiration							▪	
Titration						▪		

## 14.6 Acceleration Sensor

	Unit	Value
Type: LIS331DLH		
Quantity		1
Measuring range: 3 axes		

## 14.7 RFID Tag

	Unit	Value
Type: HTS 2048 RFIC IC		
Quantity		1
Frequency	kHz	125

## 14.8 Adjustment

### 14.8.1 Initial Adjustment

---

Type of adjustment by manufacturer in accordance with ISO 8655-1

---

On dispensing (Ex) of the selected volume (test volume), at 20 °C

---

Adjustment setting in factory settings, see Chapter "4.9.10 Parameters in the "Settings"/"Adjustment" Menu", page 29

---

## 14.9 Materials

---

### Housing

1.4404, aluminium

---

Plastic PBT | PA

---

### User display

Plastic PBT | PP

---

Float glass

---

### Tip ejector

Models with volume range 10 µL | 120 µL | 300 µL: polyvinylidene fluoride (PVDF)

---

Models with volume range 5,000 µL | 10,000 µL: polypropylene (PP)

---

### Tip cone holder

Models with volume range 10,000 µL: polyetherimide (PEI)

---

For all other models: polyamide (PA)

---

### Tip cone

Models with volume range 10 µL: polyvinylidene fluoride (PVDF)

---

Models with volume range 200 µL | 300 µL: polyetherimide (PEI)

---

Models with volume range 1,000 µL | 5,000 µL | 10,000 µL: polyphenylene sulphide (PPS)

---

### Piston

Models with volume range 3 µL | 10 µL | 20 µL: stainless steel (SS)

---

Models with volume range 100 µL (1 channel): polyphenylene sulphide (PPS)

---

Models with volume range 100 µL (multi-channel model): polyetherimide (PEI)

---

Models with volume range 200 µL | 300 µL | 1,000 µL | 5,000 µL: polyphenylene sulphide (PPS)

---

### Piston seal

Models with volume range 3 µL | 10 µL | 20 µL | 10,000 µL: fluoroelastomer (FKM)

---

Models with volume range 100 µL | 200 µL | 1,000 µL | 5,000 µL: ethylene propylene diene monomer rubber (EPDM)

---

Spring

Models with volume range 10 µL: Stainless steel (SS)

Information on interactions of materials with organic and inorganic solutions, solvents and corrosive chemicals can be provided on request.

## 14.10 Cleaning and Sterilization

### 14.10.1 Cleaning Agents, Decontamination Agents and Cleaning Methods

Approved cleaning agents and decontamination agents

Mild cleaning agent (e.g. 70% ethanol solution | 65% isopropanol solution)

Disinfectant and decontamination solution (e.g. 70% ethanol, 65% isopropanol)

For other cleaning and decontamination agents contact Sartorius.

Approved cleaning methods

Wiping the device surfaces with a slightly damp cleaning cloth

Wiping the device surfaces dry

Replacing the Safe-Cone Filters with the tweezers

### 14.10.2 Approved Sterilization Methods for Lower Assembly

	Unit	Value
Sterilization methods		
UV radiation		
Wiping with an approved decontamination agent (see Chapter 14.10.1, page 68)		
Autoclaving, only for suitable devices and only under autoclaving conditions		
Other sterilization methods or application-specific sterilization methods are possible on request		
Suitable devices for autoclaving		
1-channel model: All models		
Multi-channel model: Models with maximum volume of 10 µL   120 µL   300 µL, marked with autoclaving symbol		
Autoclaving conditions		
Autoclaving temperature, maximum	°C	121
Approved overpressure	bar	1
Duration, maximum	Minutes	20
Autoclaving in autoclaving bag		

## 14.11 Battery

	Unit	Value
Li-polymer battery with protective circuit, rechargeable		
Service life at room temperature, expected	Years	2
Capacity	mAh	350
Charging time	h	1

## 14.12 Routine test

### 14.12.1 Conditions for Aspirating the Test Liquid

Test volume ( $\mu\text{L}$ )	Immersion depth of the pipette tip during aspiration (mm)	Wait time (s)
$\leq 1$	1 – 2	1
> 1 to 100	2 – 3	1
> 100 to 1 000	2 – 4	1
> 1 000 to 20 000	3 – 6	3

### 14.12.2 Correction Factor Z ( $\mu\text{L} / \text{mg}$ ) for Routine Test

Temperature	Air pressure			
	95	100	101.3	105
Unit	Unit	Unit	Unit	Unit
$^{\circ}\text{C}$	kPa	kPa	kPa	kPa
Value	Value	Value	Value	Value
20.0	1.0028	1.0028	1.0029	1.0029
20.5	1.0029	1.0029	1.0030	1.0030
21.0	1.0030	1.0031	1.0031	1.0031
21.5	1.0031	1.0032	1.0032	1.0032
22.0	1.0032	1.0033	1.0033	1.0033
22.5	1.0033	1.0034	1.0034	1.0034
23.0	1.0034	1.0035	1.0035	1.0036
23.5	1.0036	1.0036	1.0036	1.0037

The values in the table represent the most common values. For the full table, or for formula to calculate correction factor the standard ISO 8655-6 must be considered.

## 14.13 Performance Specification

### 14.13.1 Single-Channel Models

Model	Colour of the operating button	Volume range	Maximum permissible error limits, according to ISO 8655				
			Mode**   Test volume	Systematic error*		Random error*	
				Unit	Unit	Unit	Unit
				µL	µL	%	µL
Value	Value	Value	Value	Value	Value		
LH-747021	■	0.5 – 10	P   10	1.0	0.100	0.4	0.040
			P   5	1.2	0.060	0.7	0.035
			P   1	3.0	0.030	2.0	0.020
			P   0.5	8.0	0.040	4.0	0.020
			D   1	6.0	0.060	7.0	0.070
LH-747041	■	5 – 120	P   120	0.5	0.60	0.15	0.18
			P   60	0.7	0.42	0.23	0.14
			P   12	2.0	0.24	1.0	0.12
			P   5	5.5	0.275	2.5	0.125
			D   12	4.0	0.48	4.0	0.48
LH-747061	■	10 – 300	P   300	0.6	1.80	0.15	0.45
			P   150	0.6	0.90	0.2	0.30
			P   30	2.0	0.60	0.8	0.24
			P   10	6.0	0.60	2.4	0.24
			D   30	3.0	0.90	3.0	0.90

\* The listed values for the systematic and random errors are valid under the following conditions:

- Use of the listed models with suitable Sartorius Optifit non-sterile tips
- Determination of the values under strictly controlled conditions during type examinations in accordance with ISO 8655

The values for the systematic and random errors listed in ISO 8655 are valid under the following conditions:

- Use of the listed models with other Sartorius tips

\*\* P = Pipetting mode | D = mode for Multi Dispensing

Model	Colour of the operating button	Volume range	Maximum permissible error limits, according to ISO 8655				
			Mode**   Test volume	Systematic error*		Random error*	
				Unit	Unit	Unit	Unit
			µL	µL	%	µL	%
Value	Value	Value	Value	Value	Value		
LH-747081	■	50 - 1000	P   1000	0.45	4.5	0.15	1.5
			P   500	0.6	3.0	0.2	1.0
			P   100	2.0	2.0	0.5	0.5
			P   50	4.0	2.0	1.0	0.5
			D   100	2.5	2.5	2.0	2.0
LH-747101	■	100 - 5000	P   5000	0.5	25	0.15	7.5
			P   2500	0.7	17.5	0.2	5
			P   500	1.6	8	0.4	2
			P   100	8.0	8	2.0	2
			D   500	2.4	12	2.4	12
LH-747111	■	500 - 10000	P   10000	0.6	60	0.2	20
			P   5000	0.9	45	0.3	15
			P   1000	3.0	30	0.6	6
			P   500	7.0	35	1.2	6
			D   1000	4.0	40	2.4	24

\* The listed values for the systematic and random errors are valid under the following conditions:

- Use of the listed models with suitable Sartorius Optifit non-sterile tips
- Determination of the values under strictly controlled conditions during type examinations in accordance with ISO 8655

The values for the systematic and random errors listed in ISO 8655 are valid under the following conditions:

- Use of the listed models with other Sartorius tips

\*\* P = Pipetting mode | D = mode for Multi Dispensing

## 14.13.2 Multi-Channel Models

Model	Number of Channels	Colour of the operating button	Volume range	Maximum permissible error limits, according to ISO 8655					
				Mode**   Test volume		Systematic error*		Random error*	
				Unit	Unit	Unit	Unit	Unit	Unit
				µL	µL	%	µL	%	µL
Value	Value	Value	Value	Value	Value				
LH-747321   LH-747421	8   12	■	0.5 – 10	P   10		1.2	0.120	0.5	0.050
				P   5		1.5	0.075	0.8	0.040
				P   1		4.0	0.040	3.0	0.030
				P   0,5		10.0	0.050	6.0	0.030
				D   1		12.0	0.120	15.0	0.150
LH-747341   LH-747441	8   12	■	5 – 120	P   120		0.6	0.72	0.3	0.36
				P   60		0.8	0.48	0.4	0.24
				P   12		2.5	0.30	1.67	0.20
				P   5		6.0	0.30	4.0	0.20
				D   12		4.5	0.54	8.0	0.96
LH-747361   LH-747461	8   12	■	10 – 300	P   300		0.6	1.80	0.2	0.60
				P   150		0.8	1.20	0.3	0.45
				P   30		2.33	0.70	1.0	0.30
				P   10		8.0	0.80	3.0	0.30
				D   30		3.33	1.00	6.0	1.80

\* The listed values for the systematic and random errors are valid under the following conditions:

- Use of the listed models with suitable Sartorius Optifit non-sterile tips
- Determination of the values under strictly controlled conditions during type examinations in accordance with ISO 8655

The values for the systematic and random errors listed in ISO 8655 are valid under the following conditions:

- Use of the listed models with other Sartorius tips

\*\* P = Pipetting mode | D = mode for Multi Dispensing



Model	Number of Channels	Colour of the operating button	Volume range	Maximum permissible error limits, according to ISO 8655					
				Mode**   Test volume		Systematic error*		Random error*	
				Unit	Unit	Unit	Unit	Unit	Unit
				µL	µL	%	µL	%	µL
Value	Value	Value	Value	Value	Value				
LH-747391   LH-747491	8   12	■	50 - 1200	P   1200	0.6	7.2	0.2	2.4	
				P   600	1.0	6.0	0.3	1.8	
				P   120	2.5	3.0	1.0	1.2	
				P   50	8.0	4.0	2.0	1.0	
				D   120	3.33	4.0	3.33	4.0	

\* The listed values for the systematic and random errors are valid under the following conditions:  
 – Use of the listed models with suitable Sartorius Optifit non-sterile tips  
 – Determination of the values under strictly controlled conditions during type examinations in accordance with ISO 8655

The values for the systematic and random errors listed in ISO 8655 are valid under the following conditions:  
 – Use of the listed models with other Sartorius tips

\*\* P = Pipetting mode | D = mode for Multi Dispensing

## 14.14 Speed Table

### 14.14.1 1-channel Pipette

		10 µL	120 µL	300 µL	1,000 µL	5,000 µL	10,000 µL
Speed*	Unit	Value	Value	Value	Value	Value	Value
1	s	2.5	6.0	7.7	10.1	10.2	10.2
2	s	1.8	4.2	5.3	7.4	7.4	7.4
3	s	1.3	2.9	3.7	5.4	5.4	5.4
4	s	1.0	2.1	2.7	3.8	3.8	3.8
5	s	0.8	1.5	1.9	2.8	2.7	2.9
6	s	0.6	1.1	1.4	1.9	1.8	2.2
7	s	0.5	0.9	1.1	1.2	1.1	1.7
8	s	0.4	0.7	0.9	0.8	0.8	1.3
9	s	0.3	0.6	0.8	0.6	0.6	0.9

\* The speed is measured in pipetting mode with maximum volume. The speed is graduated from 1 (slow) to 9 (fast). In all main operating modes, the speed can be set separately for aspiration and dispensing.

### 14.14.2 Multi-channel Pipette

		10 µL	120 µL	300 µL	1,200 µL
Speed*	Unit	Value	Value	Value	Value
1	s	2.5	6.1	5.4	6.1
2	s	1.8	4.4	3.9	4.4
3	s	1.3	3.3	2.9	3.3
4	s	1.0	2.4	2.1	2.5
5	s	0.8	1.8	1.6	1.9
6	s	0.6	1.4	1.2	1.4
7	s	0.5	1.1	1.0	1.1
8	s	0.4	0.9	0.8	0.9
9	s	0.3	0.7	0.7	0.7

\* The speed is measured in pipetting mode with maximum volume. The speed is graduated from 1 (slow) to 9 (fast). In all main operating modes, the speed can be set separately for aspiration and dispensing.

# 15 Accessories and Consumables

## 15.1 Accessories

This table contains an excerpt of the accessories that can be ordered. For information on further items, contact Sartorius.

Item	Quantity	Order number
Charging devices		
Charging stand for 1 pipette with universal charger	1	730981
Charging carousel for 4 pipettes with universal charger	1	730991
USB charging device	1	LH-735001
Storage option (non-charging)		
Pipette stand for all Sartorius pipettes	1	725620
Holder for 1 pipette	1	LH-727640
Reagent container (capacity 120 mL)	1	783500

## 15.2 Consumables

This table contains an excerpt of the consumables that can be ordered. For information on further items, contact Sartorius.

Item	Quantity	Order number
Optifit Tips, 0.1-10 µL, single rack	1	790010
Optifit Tips, 0.1-10 µL, refill pack	1	790013
Optifit Tips, 0.1-10 µL, refill tower	1	790011
Safetyspace® Filter-Tip, 0.1-10 µL, single rack	1	790011F
Safetyspace® Filter-Tip, 0.2-120 µL, single rack	1	790101F
Safetyspace® Filter-Tip, 50-1,000 µL, single rack	1	791001F
Extended Standard Tip, 0.1-10 µL, single rack	1	783210
Extended Standard Tip, 50-1,200 µL, single rack	1	791210
Extended Filter Tips 0.1-10 µL, single rack	1	783201
Extended Filter Tips 10-1,000 µL, single rack	1	LH-XF781001
Safe-Cone Filter, standard, Ø6.73 mm	50	721005
Safe-Cone Filter, standard, Ø5.33 mm	50	721006
Safe-Cone Filter, standard, Ø3.15 mm	50	721007
Safe-Cone Filter, standard, Ø2.51 mm	50	721008
Safe-Cone Filter, standard, Ø1.83 mm	50	721014
Tweezers for filter replacement	1	721009

## 16 Sartorius Service

Sartorius Service would be happy to help with any queries about the device. For information on the Service addresses, Service services and on the local contact, see the Sartorius website ([www.sartorius.com](http://www.sartorius.com)).

If you have queries about the system or are getting in touch in the event of malfunctions, have the device information to hand and give it to Sartorius Service, e.g. serial number, hardware, firmware, configuration. To do this, note the information on the manufacturer's ID label and in the "Settings"/"Information" menu.

## 17 Conformity of the Device

In the following, the compliance of the device with the designated directives or standards is declared.

### EU/EC Declaration of Conformity

<b>Manufacturer</b>	Sartorius Biohit Liquid Handling Oy Laippatie 1, 00880 Helsinki, Finland	
	We hereby declare under our sole responsibility that the following product	
<b>Name</b>	Picus® 2	
<b>Modell(s)</b>	LH-747021, LH-747041, LH-747061, LH-747081, LH-747101, LH-747111, LH-747321, LH-747341, LH-747361, LH-747391, LH-747421, LH-747441, LH-747461, LH-747491	
<b>Device type</b>	Electronic pipette	
	in the form as delivered fulfils all the relevant provisions of the following European Directives – including any amendments valid at the time this declaration was signed – and meets the applicable requirements of the harmonized European Standards including any amendments valid at the time this declaration was signed – listed below:	
<b>2014/53/EU</b>	Directive on Radio Equipment EN 300 328 V2.2.2 EN 301 489-17 V3.2.4	
<b>2014/30/EU</b>	Directive on Electromagnetic Compatibility EN 61326-1:2013 EN 301 489-1 V1.9.2	
<b>2011/65/EU</b>	Restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS) (incl. (EU) 2015/863) EN IEC 63000:2018	
<b>Accessories</b>	Charging Stand	730981
	Charging Carousel	730991
	Picus® AC Adapter	LH-735001
	Sartorius Pipetting mobile application	

## UK Declaration of Conformity

<b>Manufacturer</b>	Sartorius Biohit Liquid Handling Oy Laippatie 1, 00880 Helsinki, Finland	
	We hereby declare under our sole responsibility that the following product	
<b>Name</b>	Picus® 2	
<b>Modell(s)</b>	LH-747021, LH-747041, LH-747061, LH-747081, LH-747101, LH-747111, LH-747321, LH-747341, LH-747361, LH-747391, LH-747421, LH-747441, LH-747461, LH-747491	
<b>Device type</b>	Electronic pipette	
	in the form as delivered fulfils all the relevant provisions of the following UK Regulations and meets the applicable requirements of the UK Designated Standards including any amendments valid at the time this declaration was signed listed below:	
	The Radio Equipment Regulations 2017 UK Statutory Instruments 2017 No. 1206 EN 300 328 v. 2.2 EN 301 489-17 V3.2.4	
	The Electromagnetic Compatibility Regulations 2016 UK Statutory Instruments 2016 No. 1091 BS EN 61326-1:2013 EN 301 489-1 V1.9.2	
	The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012 (RoHS) UK Statutory Instruments 2012 No. 3032 BS EN 50581:2012, BS EN 61010-1:2010	
<b>Accessories</b>	Charging Stand	730981
	Charging Carousel	730991
	Picus® AC Adapter	LH-735001
	Sartorius Pipetting mobile application	

## ANATEL

The device contains ANATEL approved module. The approval number is 00857-21-05903.

## IMDA Standards



The device complies with IMDA standards: DA105282.

## 17.1 Compliance Information and Information for Users for Radio Equipment

### 47 CFR 2.1077

This device complies with part 15 of the FCC Rules. Operation is subject to the following two conditions:

- (1) This device may not cause harmful interference, and
- (2) this device must accept any interference received, including interference that may cause undesired operation.”

### 47 CFR 15.21

Changes or modifications not expressly approved by the party responsible for compliance could void the user’s authority to operate the equipment.

### 47 CFR 15.105

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 is connected.
- Consult the dealer or an experienced radio/TV technician for help.”

## 18 Trademark Information

### 18.1 Bluetooth®

Bluetooth® is a registered trademark of Bluetooth SIG, Inc..

Sartorius Liquid Handling oy  
Tulppatie 1  
00880 Helsinki, Finland

Lhinfo.finland@sartorius.com  
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:

06 | 2024

© 2024

Sartorius Liquid Handling oy  
Tulppatie 1  
00880 Helsinki, Finland

NT | Publication No.: WPI6002-e240804