



Thermo Scientific CW3 Cell Washer

Instruction Manual

50144259-i • 12 / 2021

**KONFORMITÄTSERKLÄRUNG
DECLARATION OF CONFORMITY**

Name und Anschrift des Herstellers und des Bevollmächtigten
für die Zusammenstellung der relevanten technischen Unterlagen:

*Name and address of the manufacturer and of the authorized
representative to compile the relevant technical documentation:*

**Thermo Electron LED GmbH
Zweigniederlassung Osterode
Am Kalkberg
37520 Osterode am Harz
Germany**

Gegenstand der Erklärung / *Object of the declaration:*

Beschreibung /description	: Zentrifuge mit Zubehör / Centrifuge with accessories
Marke / brand	: Thermo Scientific
Modellbezeichnung / model name	: CW3
Modell Nr. / model no.	: 75007405
Gültig ab Equipmentnr. Valid from equipment no.	: O63458

mit allen einschlägigen Bestimmungen der Richtlinie über In-vitro-Diagnostika 98/79/EG in Übereinstimmung ist.
is in conformity with all relevant terms of directive for in vitro diagnostic medical devices 98/79/EC.

Die Maschine ist auch in Übereinstimmung mit den Schutzziele der Maschinenrichtlinie 2006/42/EG, der Niederspannungsrichtlinie 2014/35/EU und der Richtlinie 2014/30/EU über elektromagnetische Verträglichkeit.
The machinery is in accordance with the protection goals for the directives machinery 2006/42/EC, low voltage 2014/35/EU and electromagnetic compatibility 2014/30/EC.

Der oben beschriebene Gegenstand der Erklärung erfüllt auch die Vorschriften der Richtlinie 2011/65/EU des Europäischen Parlaments und des Rates vom 8. Juni 2011 zur Beschränkung der Verwendung bestimmter gefährlicher Stoffe in Elektro- und Elektronikgeräten. Die alleinige Verantwortung für die Ausstellung dieser Konformitätserklärung trägt der Hersteller.

The object of the declaration described above is also in conformity with Directive 2011/65/EU of the European Parliament and of the Council of 8 June 2011 on the restriction of the use of certain hazardous substances in electrical and electronic equipment.


This declaration of conformity is issued under the sole responsibility of the manufacturer.

Angewandte Normen/
Standards used:

**EN 61010-1: 2004
EN 61010-2-020: 2006
EN 61010-2-101: 2002
EN 61326-1: 2013
EN 61326-2-6: 2013**

Unterzeichnet für und im Namen von: Thermo Electron LED GmbH.
Signed for and on behalf of: Thermo Electron LED GmbH

Osterode am Harz, den 26.10.2016


Dr. Andreas Karl,
Director R&D Global Project Management

	Name	Datum	Dokument	Revision
Freigegeben	Abdullah Laaboubi	18.10.2016	50148108	01






Table of Contents

1. Preface	5
1. 1. Items Supplied	5
1. 2. Intended Use	7
1. 3. Precautions	7
1. 4. Symbols Used on the Cell Washer	10
1. 5. Symbols Used in the Manual	11
2. Technical Specifications	12
2. 1. Directives and Standards	13
2. 2. Mains Supply	13
2. 3. Location and Function of Parts	14
2. 4. Control Panel	16
2. 5. Condition-setting Panel	17
3. Transport and Set Up	19
3. 1. Before Setting Up	19
3. 2. Location	19
3. 3. Transport	19
3. 4. Level	20
3. 5. Check Mains Connection	20
3. 6. Set Up	20
3. 7. Storage	32
3. 8. Shipping	32
4. Operation	33
4. 1. Preparation	33
4. 1. 1. Injection Volume Adjustment and Pump Priming	34
4. 1. 2. Preparation of the Sample	36
4. 2. Operation Modes	37
4. 2. 1. Automatic Mode	37

4. 2. 2. Manual Mode	37
4. 2. 3. Agitation-Centrifugation.	38
4. 2. 4. Operational Sequence for Antiglobulin Test.	38
5. Maintenance	39
5. 1. Cleaning	39
5. 1. 1. Pump, Tank and Tubing.	40
5. 1. 2. Rotor, Distributor and Bowl	40
5. 1. 3. Chamber, Splash Guard, Drain Cover and Door Stopper Parts	41
5. 2. Preventive Maintenance	42
5. 3. Shipping and Disposal	42
6. Troubleshooting	43
6. 1. Mechanical Emergency Door Release	43
6. 2. Error Codes	44
6. 3. Troubleshooting when no Error Code is indicated	45
Decontamination Information Certificate	46
Chemical Compatibility	47
Index	50



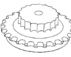

1. Preface

1. 1. Items Supplied

Article No.	Description	Quantity	
	Thermo Scientific CW3 cell washer	1	
S402776A	Bowl assembly	1	
S413259C	Tank (5 L)	1	
4744346	10 mm dia. tube	3 m	
75000015	Tube connector	1	
S4011034	18 mm dia. tube	2.5 m	
480270	Test tubes (12 mm in diameter and 75 mm in length)	50 (1 box)	
483719	Grease	1	
8046004	Hex wrench	1	
S411107	Motor guard plate	1	
4666354	<u>Hose band</u> Small	1	
4666357	Medium	1	
4666355	Large	1	
S413181A	L-shaped adapter	1	
480879	D10 adapter	25	
	Power cord	1	
	Instruction manual	1	
	Storage medium with instruction manuals	1	
	Quick installation guide	1	

Rotor Options			
75000020	24 place rotor assembly with distributor		
75000021	12 place rotor assembly with distributor		

Optional Accessories			
75000026	1 000 ml glass bottle with sensor		
75007410	Modified saline sensor kit		

Article No.	Description	Quantity	
Replacement Parts			
75000022	<u>Rotor Assembly</u> 24 places		
75000023	12 places		
75000024	<u>Distributor Assembly</u> 24 places		
75000025	12 places		

1. 2. Intended Use

The Thermo Scientific CW3 cell washer is designed to perform cell washing in multiple washing cycles using saline solution. The cell washer provides blood cells after sample separation, which can be used for further blood testing such as Antiglobulin test, ABO compatibility, Rh testing, Cross-matching and Antibody screening.

The centrifuge should always be operated by a trained individual such as a clinical laboratory technologist or a person with a similar education.

1. 3. Precautions

Carefully read the following safety precautions for a thorough understanding.

Protection is impaired if the cell washer is not used as specified in the intended use.

Follow the instructions and procedures described in this instruction manual to operate this cell washer safely.

Be sure to observe the precautions in this instruction manual and on your equipment.

It is the general obligation of the operator to make sure that the proper protective clothing is used. Mind the „Laboratory Biosafety Manual“ of the World Health Organization (WHO) and the regulations in your country.

Safety messages are indicated as follows. They are stated with signal words combined with the safety alert symbol to call your attention to items or operations that could be hazardous to you or other persons using this equipment. The definitions of signal words are as follows:

- **WARNING:** Warning notes indicate any condition or practice, which if not strictly observed, could result in serious injury or possible death.
- **CAUTION:** Caution notes indicate any condition or practice, which if not strictly observed or stopped, could result in personal injury, damage or destruction of the equipment.
- **NOTICE:** Notes indicate an area or subject of special merit, emphasizing either the product's capabilities or common errors in operation or maintenance.

Do not operate this cell washer in any manner not described in this Instruction Manual. Should you have any troubles with this cell washer, call a Thermo Fisher Scientific authorized sales/service representative.

The precautions described in this Instruction Manual are carefully developed to cover all possible risks. It is important that you are alert to unexpected incidents. Be careful when operating this cell washer.



WARNING

1. The cell washer is neither inert nor protected against explosion. Never use the cell washer in an explosion-prone environment.
2. Do not set up the cell washer in or near places where flammable gases are generated or chemicals are stored.
3. If centrifuging any hazardous materials mind the „Laboratory Biosafety Manual“ of the World Health Organization (WHO). Look on the internet page of the World Health Organization (www.who.int) for the „Laboratory Biosafety Manual“.
For materials in a higher risk group, extra safety measures have to be taken.
4. Take all necessary safety measures before using samples that are toxic or radioactive, or blood samples that are pathogenic or infectious. You use such samples at your own responsibility.
 - a. If the cell washer, rotor, or an accessory is contaminated by samples that are toxic or radioactive, or blood samples that are pathogenic or infectious, be sure to decontaminate the item according to good laboratory procedures and methods.
 - b. If there is a possibility that the cell washer, rotor, or an accessory is contaminated by samples that may impair human health (for example, samples that are toxic or radioactive, or blood samples that are pathogenic or infectious), it is your responsibility to sterilize or decontaminate the cell washer, rotor, or the accessory properly before requesting repairs from a Thermo Fisher Scientific authorized sales/service representative.
 - c. It is your responsibility to sterilize and/or decontaminate the cell washer, rotor, or parts properly before returning them to a Thermo Fisher Scientific authorized sales/service representative.
5. Do not centrifuge explosive or flammable materials or substances that could react violently.
6. Because liquid samples, saline, and other substances are used in this cell washer, your cell washer must be grounded properly.
7. To avoid electrical shocks, do not handle the power cord or turn on or off the POWER switch with wet hands.
8. The cell washer itself may move if the rotor fails during high-speed rotation. Make sure there is a 30 cm area around the cell washer that will allow for such movement and do not allow individuals to enter into that area during operation. Also do not place hazardous objects such as flammable or explosive materials on top of the cell washer or in the surrounding area.
9. While the rotor spins, never unlock the door.
10. Repairs, disassembly, and other modifications to the cell washer are strictly prohibited unless performed by a Thermo Fisher Scientific authorized sales/service representative.



CAUTION

1. Turn off the cell washer at the main switch. The mains plug must be freely accessible at all times.
Press STOP to shut the cell washer down.
Pull out the power supply plug or disconnect the power supply in an emergency.
2. Use the cell washer only with the rotor and distributor assembly correctly installed, and the bowl, splash guard and drain cover correctly inserted.
3. Do not move or relocate the cell washer while the rotor is spinning.
4. Do not lean on the cell washer.
5. Do not pour any liquids such as water, detergent, or disinfectant directly into the rotor chamber. If you do so, the bearings of the drive unit may corrode or deteriorate.
6. Connect the saline tank to the pump inlet connector of the cell washer and secure them correctly using the hose bands (standard accessories). Otherwise liquid may leak from the tube and penetrate into the cell washer. If you suspect that liquid is inside of the cell washer, contact a Thermo Fisher Scientific authorized sales/service representative to clean and dry the cell washer.
7. Before you operate the cell washer, remove any dropped objects and tube fragments from inside the rotor chamber.
8. Always check for corrosion and damage on the rotor before using it. Do not use a corroded or damaged rotor.
9. Be sure to use the specified test tubes with a wall thickness of at least 1 mm. When using thin test tubes such as disposable test tubes, do not use them repeatedly. If any scratches, cracks, inside warps, or other irregularities are found on the test tubes, do not use them because they cannot bear centrifugal force.
10. Be sure to load test tubes to all the holders.
11. Use one or two drops (about 50 µl) of 3-5% erythrocyte suspension as sample volume for blood cell washing. When using precipitated erythrocyte layer, use one or two drops (about 50 µl or less) per test tube. Use 80% or less of the test tube capacity as sample volume for centrifugation.
12. If you observe some abnormality in this product, stop using it immediately and contact a Thermo Fisher Scientific authorized sales/service representative. Notify the service representative of the alarm code, if displayed.
13. If the cell washer will not be used for a long time, remove the power cord from the socket.
14. For connection to a different power outlet, the power cord may need to be replaced.
15. Follow local electrical codes.
16. Depending on the magnitude, an earthquake may damage the cell washer.
If you observe any abnormality, contact a Thermo Fisher Scientific authorized sales/service representative.

NOTICE

Usually the control panel and the surface of the cell washer get warm during operation.

1. 4. Symbols Used on the Cell Washer



This symbol refers to general hazards.

CAUTION means that material damage may occur.

WARNING means that injuries or material damage or contamination may occur.

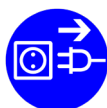


This symbol refers to biological hazards.

Observe the information contained in the instruction manual to keep yourself and your environment safe.



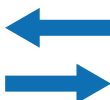
This symbol refers to information on general hazards, described within the manual.



This symbol instructs you to disconnect AC mains power before transporting or servicing the cell washer.



This symbol instructs you not to pour the open cell washer with water.



This symbol shows the inflow and draining directions of the cell washer. The arrow pointing towards the cell washer shows the inflow direction. The arrow pointing away from the cell washer shows the draining direction.



This symbol instructs you to make sure that the drain cover is installed in the cell washer door before you start the CW3 cell washer. Failure to install the drain cover raises the risk of a biological hazard when contaminated samples are used. See "5. 1. 3. 1. Drain Cover" on page 41 for removal and installation.



Caution: Federal law restricts this device to sale by or on the order of a qualified clinical facility manager or equivalent.



The CE mark states that this product meets all requirements for the European Economic Area.



This CSA mark states that this product meets all requirements for Canada and USA.



This symbol marks the manufacturer of this product.



This symbol marks the manufacturing date of this product.



This symbol marks the catalog number of this product.



This symbol marks this product as intended for use as an in vitro diagnostic medical device.

1. 5. Symbols Used in the Manual



This symbol refers to general hazards.
CAUTION means that material damage may occur.
WARNING means that injuries or material damage or contamination may occur.



This symbol refers to biological hazards.
Observe the information contained in the instruction manual to keep yourself and your environment safe.

2. Technical Specifications

Environmental Conditions	Altitudes of up to 2 000 m above sea level Max. relative humidity 80% up to 31 °C; decreasing linearly up to 50% relative humidity at 40 °C
Environmental Conditions during Storage and Shipping	Temperature: -10 °C to +55 °C Humidity: 15% to 85%
Permissible Ambient Temperature during Operation	+2 °C to +35 °C

Average Heat Dissipation	54 Wh / 184 Btu/h / 194 kJ/h
Overvoltage Category	II
Pollution Degree	2
IP	20
Running Time	Automatic: 1 – 99 sec / manual: 1 – 999 sec
Maximum Speed ² n_{max}	Centrifuge: 3 000 rpm Wash: 1 200 rpm Decant: 350-500 rpm
Maximum RCF-Value at n_{max}	Centrifuge: 1 180 x g Wash: 190 x g
Noise Level at Maximum Speed ¹	< 53 dB (A)
Maximum Kinetic Energy	0.46 kJ

Dimensions	
Height (open door / closed door)	630 mm / 410 mm
Width	370 mm
Depth	450 mm
Weight	28 kg

¹ Front Side Measurement, 1 m in front of the instrument at 1.6 m height. 3 washing-cycles each with about 35 s centrifugation.

² Mode manual centrifuge

² Mode manual decant

2. 1. Directives and Standards

Region	Directive	Standard
Europe 220-230 V, 50/60 Hz	98/79/EC In Vitro Diagnostic Medical Devices (IVD) 2014/35/EU Low Voltage (protective goals) 2006/42/EC Machinery (protective goals) 2014/30/EC Electromagnetic Compatibility (EMC) (protective goals) 2011/65/EC RoHS – Directive on the Restriction of the use of certain Hazardous Substances in electrical and electronic equipment	EN 61010-1 EN 61010-2-020 IEC 61010-2-101 EN 61326-1 Class B EN ISO 13485
USA & Canada 120 V, 60 Hz	FDA – 510(k) cleared Regulation Name: Automated Cell-Washing Centrifuge for Immunohematology Regulatory Class: 2 Product Code: KSN	CAN/CSA-C22.2 No. 61010-1-04 UL Std. No. 61010-1 CAN/CSA-C22.2 No. 61010-2-020-09- Part 2-020 IEC 61010-2-020 IEC 61010-2-101
China 220-230 V, 50/60 Hz		IEC 61010-1 IEC 61010-2-020 IEC 61010-2-101 EN 61326-1 Class B

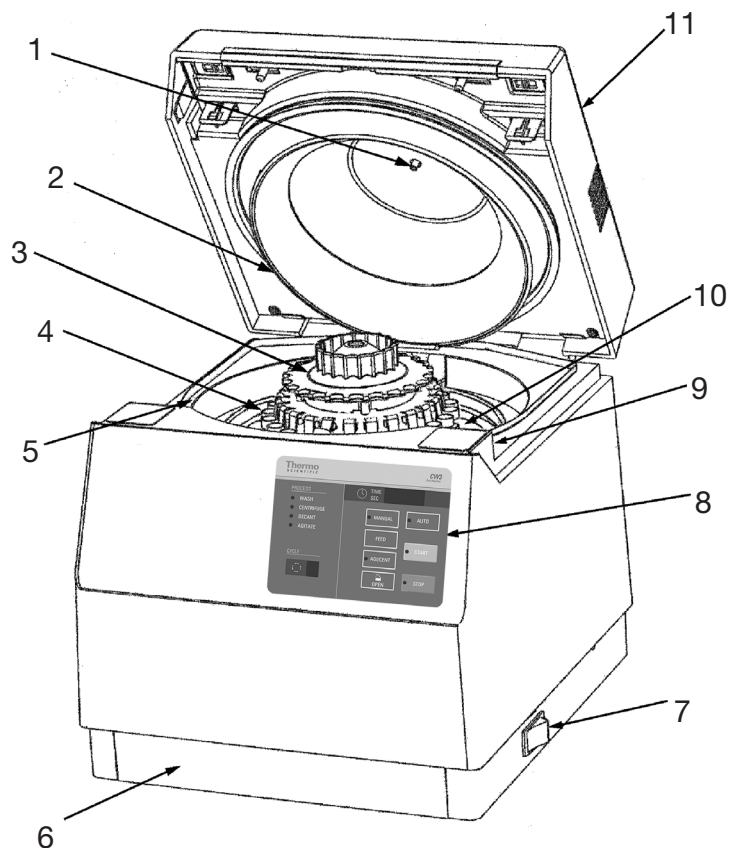
2. 2. Mains Supply

The following table contains an overview of specifications for the electrical power source. This data is to be taken into consideration when selecting the mains connection socket.

Unit	Thermo Scientific CW3 cell washer	
Article No.	75007404	75007405
Mains Voltage	120 V \pm 10%	220-30 V \pm 10%
Frequency	60 Hz	50 / 60 Hz
Rated Current	2.7 A	1.5 A
Power Consumption	135 W	135 W
Equipment Fuse	10 AT	10 AT
Building Fuse	15 AT	16 AT

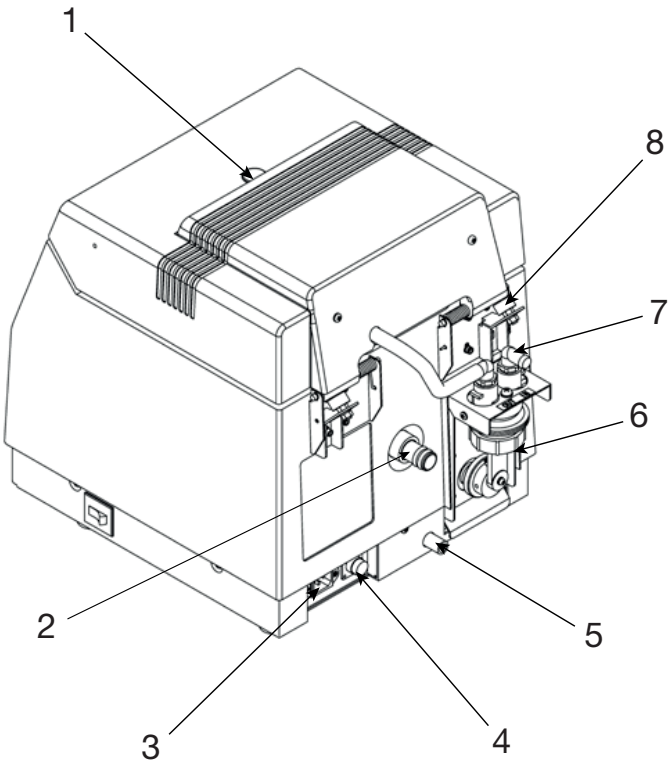
2. 3. Location and Function of Parts

Front view



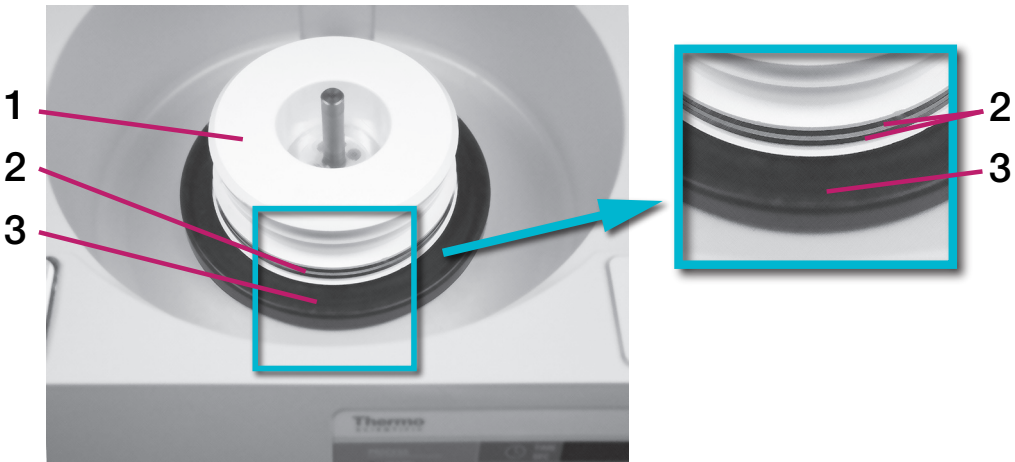
No.	Part	Description
1	Nozzle	Through this nozzle, saline is supplied from the pump to the distributor.
2	Drain cover	Receives the waste liquid decanted from the rotor and discharges it to the outside.
3	Distributor	Distributes the saline supplied from the pump to each test tube held on the rotor.
4	Rotor	A 12 or 24-sample rotor.
5	Splash guard	Designed to keep saline, waste liquid, etc. from leaking to the internal mechanism of the cell washer. Removable.
6	Condition-setting panel	See "2. 5. Condition-setting Panel" on page 17.
7	POWER switch	Switches on and off the cell washer power. It doubles as a circuit protector. It automatically switches off cell washer power when the device draws excessive currents.
8	Control panel	See "2. 4. Control Panel" on page 16.
9	Door lock lever	Keeps the door locked while the rotor is running.
10	Bowl	Rotates together with the rotor. It determines the angle at which the test-tube holders of the rotor swing out.
11	Door	Test tubes may be installed on the rotor after opening the door. The drain cover and nozzle are mounted on the bottom side of the door.

Rear view



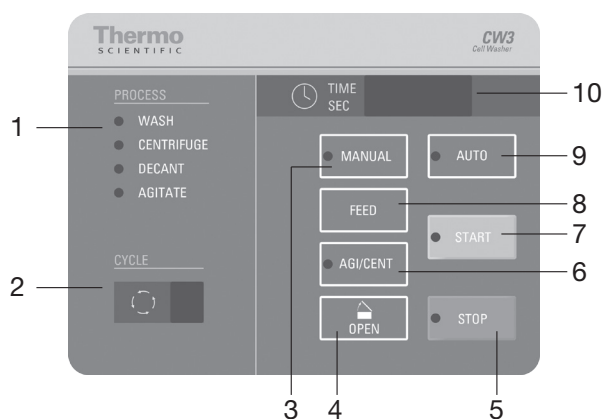
No.	Part	Description
1	View port	You can measure the rotor speed through the view port by an optical reflection type rotation meter.
2	Drain connector	Overflowing waste liquid is drained through this connector.
3	Power connector	To be connected to a power supply.
4	Sensor connector	To be connected to the sensor that detects the level of saline in the tank.
5	Auxiliary 18 mm dia. tube	The chamber is drained through this tube if the drain connector is blocked.
6	Pump	Supplies saline to the cell washer.
7	Pump inlet connector	To be connected to the saline tank by tubing.
8	Door stopper parts	Designed to limit the opening angle (60°) of the door to prevent the door opening too wide.

Top view



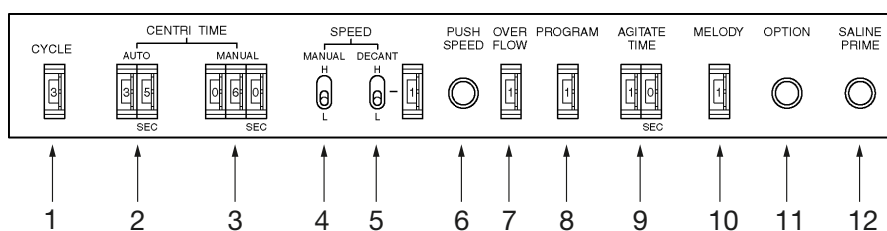
No.	Part
1	Decantation coil
2	O-ring (2x)
3	Rubber gasket

2. 4. Control Panel



No.	Name	Function
1	PROCESS LED	Shows the current process step in blood cell washing. Indicates the process setting in manual mode. While the cell washer operates, the LED of the active process step blinks.
2	CYCLE indicator	Indicates the number of washing cycles. While the cell washer operates, the CYCLE indicator shows the number of the remaining cycles.
3	MANUAL pushbutton	To manually start a single process (WASH, CENTRIFUGE, DECANT or AGITATE). The LED lights when manual mode is set.
4	OPEN pushbutton	Opens the door. If this pushbutton is pressed continuously, the cell washer does not accept any input from any of the other pushbuttons for 4 seconds (alarm buzzer sounds).
5	STOP pushbutton	Stops a process. It is also used for stopping the end buzzer and alarm buzzer. When an operation is stopped in automatic mode or during agitation/centrifugation (see No. 9), the operation can be restarted at the stopping point by pressing the START pushbutton. When the operation is stopped in manual mode, the operation is not restarted at the stopping point by pressing the START pushbutton. The STOP pushbutton LED lights when the operation is stopped. When the operation is stopped in automatic mode, this LED blinks.
6	AGI/CENT pushbutton	<p>Sets the agitation time (agitation of liquid in test tubes) and the centrifugation time (manual mode) (see the figure below). The AGI/CENT pushbutton LED lights when AGI/CENT is set. AGI/CENT operation is started by pressing the START pushbutton.</p> <p style="text-align: center;">Manual-mode set speed</p> <p>1 Agitation set time 2 Manual-mode centrifugation time</p> <p>This cell washer does not accept the START pushbutton when the agitation set time is at 0 seconds.</p>
7	START pushbutton	Starts a series of processes or a single process.
8	FEED pushbutton	In automatic mode, the FEED pushbutton causes the value of washing cycle setting to count down. In manual mode, it is used for setting the desired parameters for the current process.
9	AUTO pushbutton	Starts automatic washing cycles. The LED lights when automatic mode is set.
10	TIME indicator	Indicates the remaining time of centrifugation (in 3 digits). When an error condition occurs, the proper error code is shown. It also shows the rotor speed ($\times 10$ rpm) when PUSH SPEED on the condition-setting panel is pressed.

2. 5. Condition-setting Panel



No.	Name	Function																			
1	CYCLE setting switch	Sets the number of repetitions for the washing cycle (9 cycles at maximum).																			
2	AUTO CENTRI TIME setting switch	Sets the centrifugation time for automatic-mode operation (99 seconds at maximum). The standard centrifugation time is 35 seconds.																			
3	MANUAL CENTRI TIME setting switch	Sets the centrifugation time for manual-mode operation (999 seconds at maximum).																			
4	MANUAL SPEED setting switch	Sets one of the following rotor speeds for manual-mode centrifugation: 3,000 rpm (High) or 1,200 rpm (Low)																			
5	DECANT SPEED setting switch	<div>Sets the rotor speed for automatic- or manual-mode decantation by choosing H or L on the H/L switch and adjusting a digit between 0 and 3. Setting any digit between 4 and 9 lets the rotor default to the same speed as if 0 is set.</div> <table><tr><th rowspan="2">Switch</th><th colspan="4">Digit</th></tr><tr><th>0</th><th>1</th><th>2</th><th>3</th></tr><tr><td>H</td><td>350 rpm</td><td>400 rpm</td><td>450 rpm</td><td>500 rpm</td></tr><tr><td>L</td><td>330 rpm</td><td>370 rpm</td><td>410 rpm</td><td>450 rpm</td></tr></table> <div>Decantation is completed in a short time. Note that the above value may not necessarily match the speed indicated by the PUSH SPEED pushbutton.</div> <div>If you increase DECANT SPEED, the amount of the remaining blood cells tends to decrease.</div> <div>If you decrease DECANT SPEED, the amount of the remaining blood cells tends to increase.</div>	Switch	Digit				0	1	2	3	H	350 rpm	400 rpm	450 rpm	500 rpm	L	330 rpm	370 rpm	410 rpm	450 rpm
Switch	Digit																				
	0	1	2	3																	
H	350 rpm	400 rpm	450 rpm	500 rpm																	
L	330 rpm	370 rpm	410 rpm	450 rpm																	
6	PUSH SPEED pushbutton	Keep the PUSH SPEED pushbutton pressed when the cell washer starts spinning. The actual spinning speed of the rotor will be shown (×10 rpm) on the TIME indicator on the control panel. Example: 1200 rpm are shown as 120 on the TIME indicator (120 × 10 = 1200).																			
7	OVERFLOW setting switch	<div>When the pump is set to overflow the cell washer will overflow the tubes in the rotor with solution from the saline tank.</div> <table><tr><th>Setting</th><th>Pumping operation for overflow</th></tr><tr><td>0</td><td>No operation</td></tr><tr><td>1</td><td>Operates only during the first cycle.</td></tr><tr><td>2</td><td>Operates only during the first and the second cycles.</td></tr><tr><td>3</td><td>Operates only during the first, second and the third cycles.</td></tr><tr><td>4</td><td>Operates only during the first, second, third and the fourth cycles.</td></tr><tr><td>5</td><td>Operates only during the first, second, third, fourth and the fifth cycles.</td></tr><tr><td>6 to 9</td><td>Selectable up to 9 in the same manner.</td></tr></table>	Setting	Pumping operation for overflow	0	No operation	1	Operates only during the first cycle.	2	Operates only during the first and the second cycles.	3	Operates only during the first, second and the third cycles.	4	Operates only during the first, second, third and the fourth cycles.	5	Operates only during the first, second, third, fourth and the fifth cycles.	6 to 9	Selectable up to 9 in the same manner.			
Setting	Pumping operation for overflow																				
0	No operation																				
1	Operates only during the first cycle.																				
2	Operates only during the first and the second cycles.																				
3	Operates only during the first, second and the third cycles.																				
4	Operates only during the first, second, third and the fourth cycles.																				
5	Operates only during the first, second, third, fourth and the fifth cycles.																				
6 to 9	Selectable up to 9 in the same manner.																				

No.	Name	Function	
8	PROGRAM switch	Sets the operation program.	
		Setting	Program (No setting more than 1)
		0	5-second centrifugation is not performed during the final process of the final cycle in automatic mode.
		1	5-second centrifugation is performed during the final process of the final cycle in automatic mode.
		2	<p>You can add conditions to the centrifugation process in manual mode:</p> <ul style="list-style-type: none">▪ “3000 rpm for 15 seconds”: “H” is displayed on the CYCLE indicator and “15” is displayed in the TIME indicator.▪ “1200 rpm for 60 seconds”: “L” is displayed on the CYCLE indicator and “60” is displayed in the TIME indicator. <p>See the following procedure:</p> <ol style="list-style-type: none">1. Press MANUAL. The condition displayed on the condition-setting panel is performed.2. Press FEED. Centrifugation (3000 rpm for 15 seconds) is performed.3. Press FEED again. Centrifugation (1200 rpm for 60 seconds) is performed. <p>5-second centrifugation is not performed during the final process of the final cycle in automatic mode.</p>
(Setting the dial to any value between 3 and 9 launches the same operation program as dial setting 0.)			
9	AGITATE TIME setting switch	Sets the agitation time (99 seconds at maximum).	
10	MELODY setting switch	Sets the end buzzer.	
		Setting	Tunes
		0	No sound
		1	Electronic beep (three beeps)
		2	Oh Susanna
		3	My Bonnie
		4	My Old Kentucky Home
		5	De Camptown Races
		6, 7, 8 or 9	Electronic beep (two beeps)
You can turn down the sound level by covering the sound outlet port with adhesive tape (this port is above the right side of the CYCLE setting switch).			
11	OPTION pushbutton	This is a pushbutton intended for saline calibration purposes.	
12	SALINE PRIME pushbutton	Use for pump priming at the time of initial operation. Press this pushbutton to operate the pump for saline priming.	
		This pushbutton is ineffective while the cell washer is in operation.	

3. Transport and Set Up

3. 1. Before Setting Up

1. Check the cell washer and the packaging for any shipping damage. Inform the shipping company and Thermo Fisher Scientific immediately if any damage is discovered.
2. Remove the packaging.
3. Check if the items supplied are complete. „Items Supplied“ on page 5.
If the items supplied are incomplete, please contact Thermo Fisher Scientific service.

3. 2. Location

WARNING The cell washer is neither inert nor protected against explosion. Never use the cell washer in an explosion-prone environment.

WARNING UV rays reduce the stability of plastics. Do not subject the cell washer, rotor and plastic accessories to direct sunlight.

The cell washer is only to be operated indoors.

The set-up location must fulfill the following requirements:

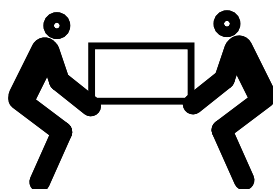
- Set up the cell washer in a room where the ambient temperature is at a constant 5 to 35 °C.
- A safety zone of at least 30 cm must be maintained around the cell washer.
Make sure that persons stay out of this safety zone while the device is centrifuging.
- The supporting structure must be stable and free of resonance.
- The supporting structure must be suitable for horizontal setup of the cell washer.
- The set-up location must be well-ventilated at all times.
- The cell washer is not to be exposed to heat and strong sunlight.

3. 3. Transport

Due to its weight, the cell washer should be carried by at least two people. Always lift the cell washer at both sides.

To prevent the risk of injuries at least two people should lift and carry the cell washer by holding it at the bottom from opposite sides.

WARNING Always lift the cell washer from both sides. Never lift the cell washer by its front panel, its back panel or at its door. Always remove the rotor before moving the cell washer.



Transport the cell washer and accessories upright within the associated packaging, if possible.

NOTICE The original cell washer packaging is a one way, disposable package. We recommend that you retain only the two large styrofoam pieces for setting up the cell washer ("3. 6. Set Up" on page 20) and properly dispose of the rest of the disposable package. Contact a shipping company or customer service for transport. Always remove the rotor before moving the cell washer. If you do not remove the rotor you may damage the cell washer drive or cell washer spindle.

3. 4. Level



CAUTION

If the cell washer is not correctly leveled, imbalance may occur and the cell washer can be damaged.
Do not place anything under the cell washer feet to level the cell washer.

Place the cell washer on horizontal and level supporting structures or benching.

Measure horizontal levelness when moving the cell washer to a new location.

3. 5. Check Mains Connection

NOTICE

Plug the cell washer into properly grounded electrical sockets only.

1. Turn off the power supply switch.
2. Check whether the power supply cable complies with the safety standards of your country.
3. Make sure that the voltage and frequency correspond to the figures on the rating plate.
4. Make sure that the power supply cable is plugged in properly.

3. 6. Set Up

NOTICE

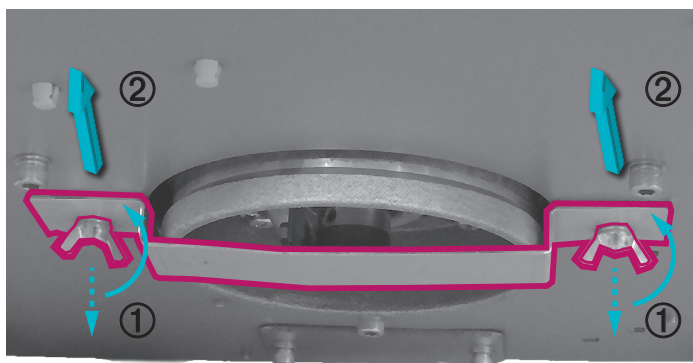
Setting up the CW3 Cell Washer is easier when done by 2 people.

1. Install motor guard plate

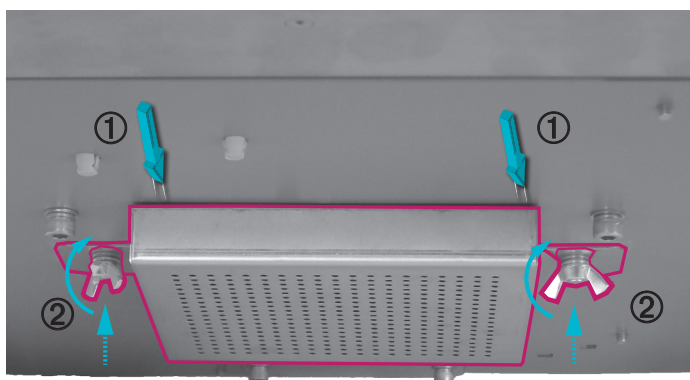
1. To remove the motor holder, put the unit on the left side (as seen from the front). Use the styrofoam blocks as support.



2. ① Loosen the wing nuts counter-clockwise.
3. ② Remove the motor holder at the bottom of the unit.



4. ① Install the motor guard plate.
5. ② Tighten the wing nuts.



6. Return the unit to the upright position.

NOTICE Keep the removed motor holder for future transportation.

2. Connect power cable

Connect the power cord to the power connector at the rear of the cell washer.

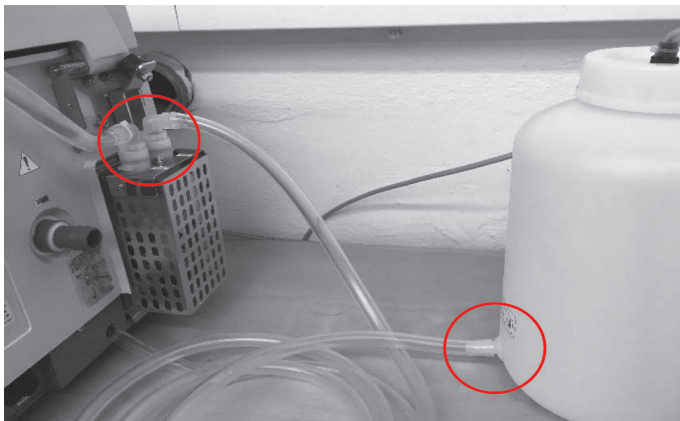
Plug the power cord in the socket.

NOTICE Plug the cell washer in grounded electrical sockets only.



3. Connect saline tank

1. Connect the bottom outlet of the saline tank to the inlet of the device using the 10 mm dia. tube. Secure both ends with the small and medium hose bands.



2. Connect the 18 mm dia. tube to the drain port in order to direct waste liquid to a tank or sink.
Optional: Use the L-shaped adapter if necessary. Secure it with the largest hose band.



3. Secure the other tubes with the remaining hose bands.



4. Connect sensor cable

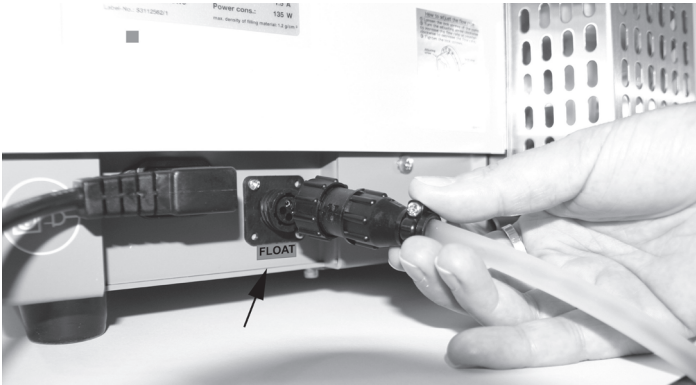
Connect the liquid sensor cable from the top of the tank to the FLOAT connector at the rear of the cell washer (see the highlighted sensor cable on the picture in „17. Connect saline source“ on page 30).

5. Switch on cell washer

Switch on the power supply switch.

The liquid level sensor in the tank activates. Error code “E2” (the saline tank needs to be refilled) is indicated and the buzzer sounds during the power-on. The buzzer stops as STOP is pressed or on its own after a short time.

6. Open door



Press OPEN and open the door.

7. Switch off cell washer

Switch off the power supply switch.

8. Remove packaging

Remove the packaging (4) from the rotor chamber.



(1) Fasteners, (2) drain cover, (3) splash guard, (4) packaging; (5) rubber gasket

9. Remove drain cover

1. Slide the two fasteners outwards to the open position.

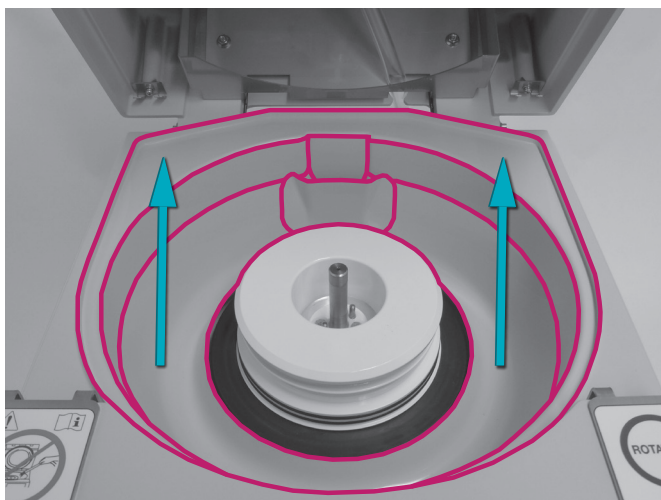


2. Pull the drain cover and fully remove it from the unit.



10. Remove splash guard

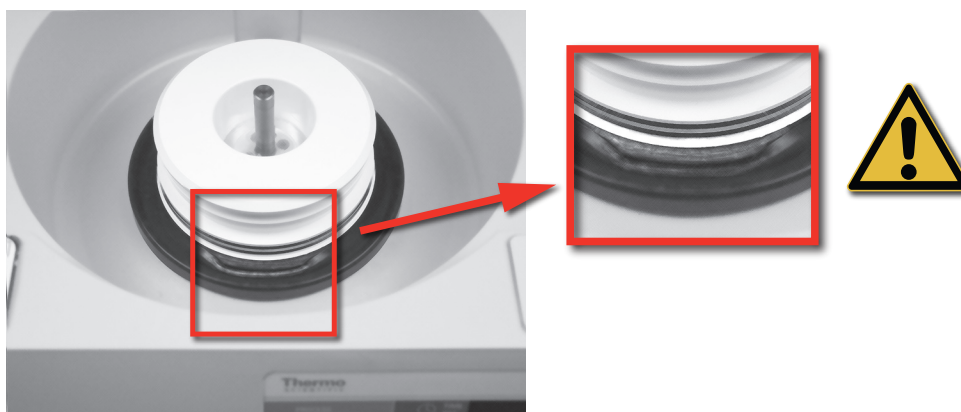
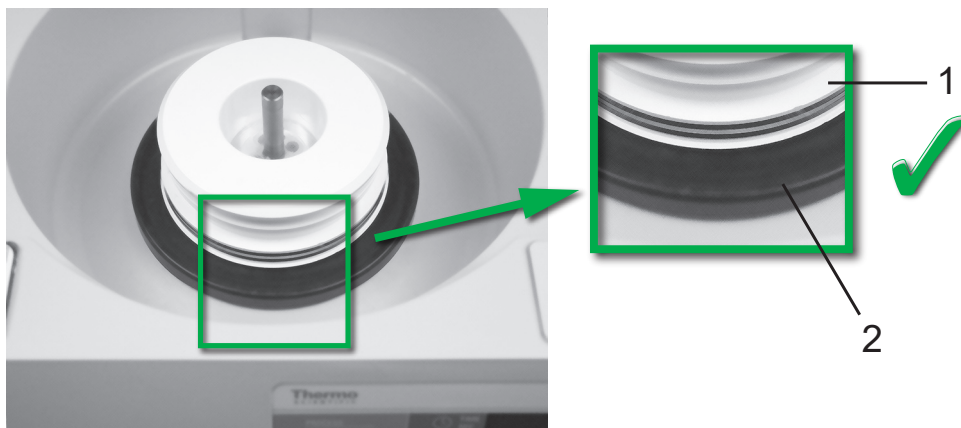
Lift and take out the splash guard.



11. Check rubber gasket

1. Check whether the white decantation coil (1) rotates freely without rubbing against the black rubber gasket (2).
2. Inspect the black rubber gasket. It should be seated without warping (as shown in the bottom picture) all around its circumference, as shown in the top picture.

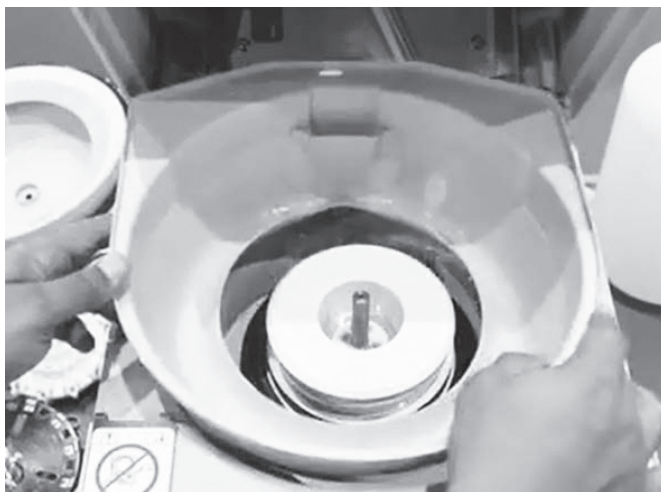
To correct the seating, push the decantation coil in one direction and the rubber gasket in the opposite direction. Continue like this until the rubber gasket is properly seated and flat against the bottom of the chamber, as shown in the top picture.



If the rubber gasket is not in position, fluids can damage the cell washer.

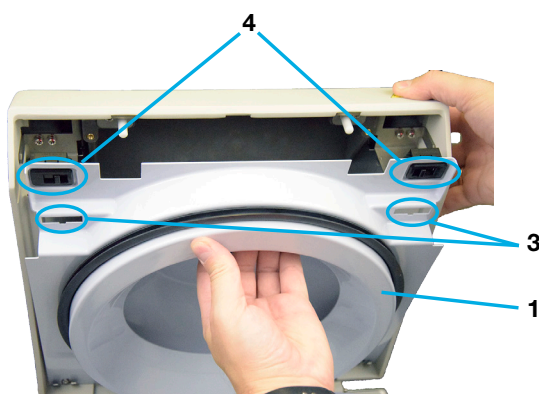
12. Install splash guard

Put the splash guard into the rotor chamber.



13. Install drain cover

1. Align the bottom end of the drain cover (1) with the hinges at the rear side of the lid.



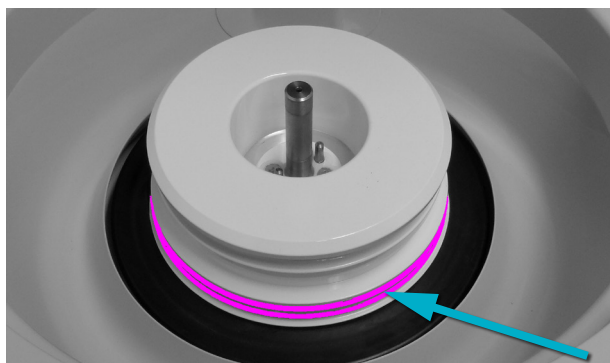
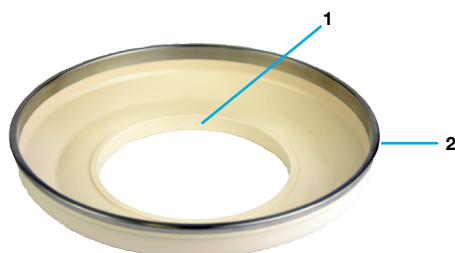
2. Align the lid hooks (2) with the slots (3) of the drain cover and push the drain cover firmly into the lid. The drain cover has to be pushed beyond the springs of the hooks. Press the top of the drain cover firmly into place until the two fasteners (4) engage with a clicking sound.



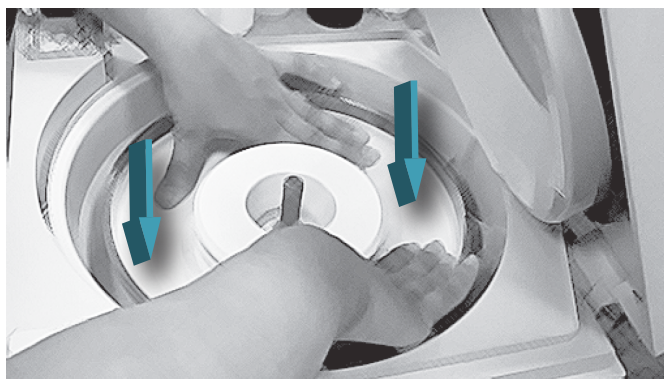
CAUTION When installing the drain cover, check that the nozzle at the center of the door is put through the hole of the drain cover. Install the drain cover correctly.

14. Install bowl

1. Apply grease (483719) delivered with the device onto the inner rim (1) of the bowl (2) and both o-rings of the decantation coil.



2. Center the bowl around the white decantation coil and push it with force into end position.



3. Push the bowl all the way down until the edge of the decantation coil becomes flush with the inner rim of the bowl. Check the position of the bowl by passing fingers over the surface. The bowl and the decantation coil must be flush with each other without any remaining step.





- ① – Bowl
- ② – Decantation coil
- ③ – O-rings

NOTICE The bowl must be installed correctly and aligned with the decantation coil before performing a test run. Make sure that the decantation coil and bowl are flush with each other by passing fingers over the seam between both parts.

If the bowl is not installed correctly, the door lock will not work, the door cannot be opened, the bowl will rub against the drain cover and the cell washer will be out of order. If you set the bowl correctly, the tube holders will touch the metal part of the bowl when the rotor is installed.

Turn the bowl by hand to verify that it rotates freely. If the bowl cannot be turned freely, the rubber gasket may not be mounted properly.



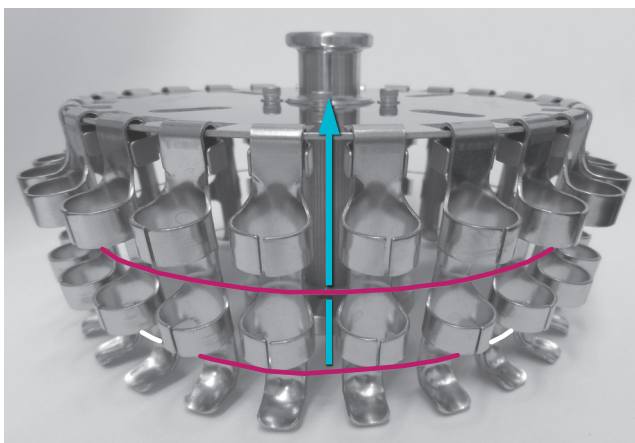
CAUTION

Error message E14 or E16 may be displayed if the bowl is not installed correctly („11. Check rubber gasket“ on page 25).

15. Install rotor

NOTICE The rotor is just set on the drive shaft. The rotor is not screwed down or otherwise tightened to the drive shaft.

1. Remove the rubber bands from the rotor.



2. Set the rotor on the drive shaft.
3. Turn the rotor about a quarter so that the coupling pins engage with the rotor shaft holes.
4. Turn the top of the rotor clockwise and counterclockwise to check that the coupling pins are engaged with the rotor shaft holes (no free turn) after setting.



CAUTION *If the rotor is not set correctly, the rotor holder may come off the bowl and overswing.*

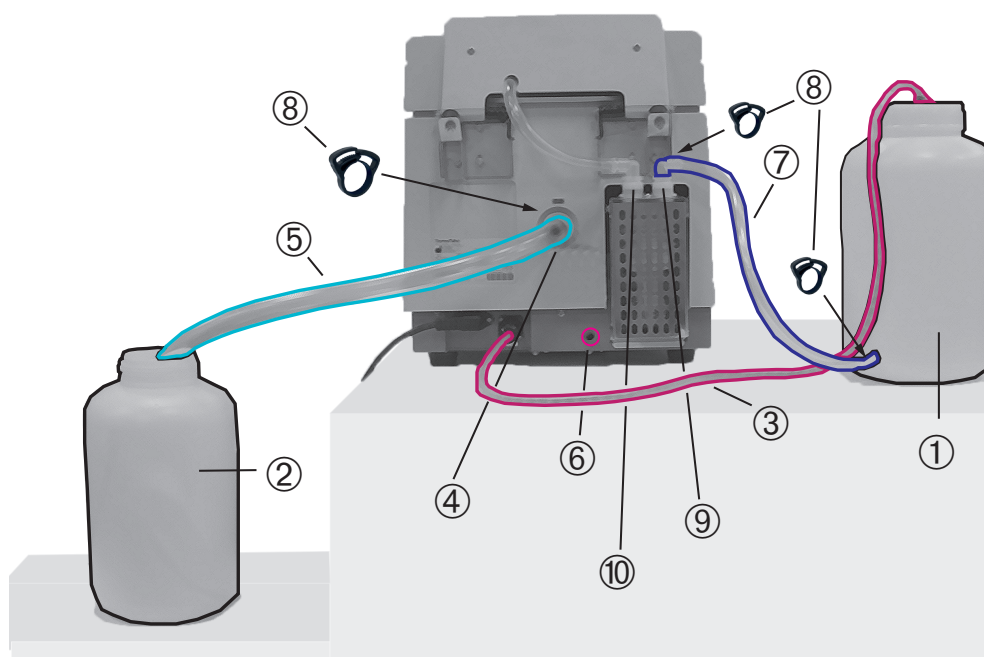
16. Install distributor

Set the distributor on the top of the rotor.



Make sure that the coupling pins at the bottom of the distributor are properly engaged with the holes at the top of the rotor.

17. Connect saline source



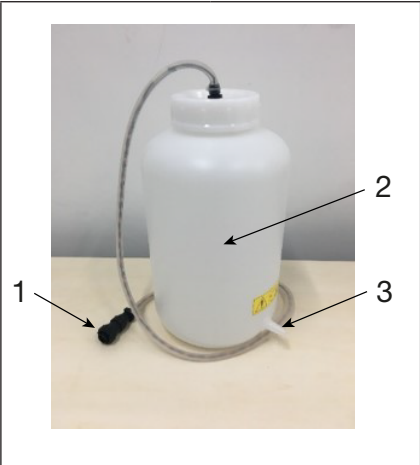
①	Saline tank
②	Waste drain or bottle (not included in items supplied)
③	Sensor cable
④	DRAIN connector
⑤	18 mm dia. tube
⑥	Gray outlet
⑦	10 mm dia. tube
⑧	Hose bands (large, medium, small)
⑨	Pump inlet connector
⑩	Pump outlet connector

CAUTION Do not set the saline tank at a place higher than the cell washer. ① The 10 mm dia. tube connected to the cell washer must be positioned higher than the other end connected to the saline tank. Otherwise the saline will overflow the cell washer and affect the effectiveness of its operation.

CAUTION If the tubing is not secured using the hose bands liquid may leak from the tube and penetrate into the cell washer. If you suspect that liquid may have leaked into the inner space of the cell washer, contact a Thermo Fisher Scientific authorized sales/service representative for cleaning and drying the cell washer.

Connect 5L saline tank to CW3 unit

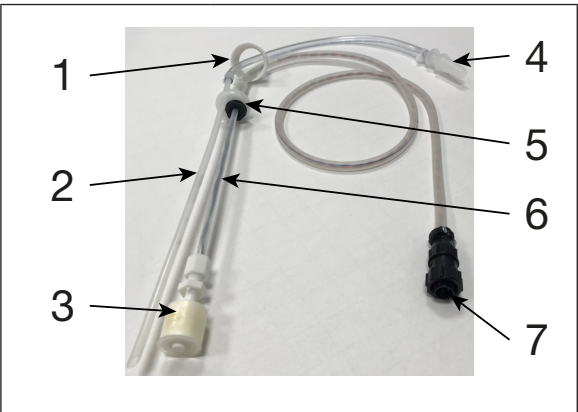
1. Place the 5L saline tank at position not higher than the cell washer unit.
2. Insert the liquid sensor into the saline tank.
3. Fill the saline tank with saline to the level higher than the liquid sensor.
4. Close the lid of saline tank.
5. Apply a thin layer of grease (483719) to the inner side of the 10mm dia. inlet tube (7) and connect it to the bottom outlet of the saline tank and the pump inlet of the cell washer (9).
Secure the tube at both ends with the hose bands.
6. Connect the sensor cable to the connector at the back of cell washer.



No.	Part
1	Sensor cable
2	Liquid sensor inside tank
3	Outlet

Connect saline cube to CW3 unit using the modified saline sensor kit (75007410)

1. Place the saline cube at position not higher than the cell washer unit, with opening facing upwards.
2. Insert the liquid sensor and inlet tube together into the saline cube.
3. Loosen the cord lock and put the liquid sensor to the bottom of the saline cube by moving up/down the sensor tube.
4. Fasten the cord lock to fix the position of liquid sensor.
5. Adjust the inlet tube to the bottom of the saline cube.
6. Tighten the screw cap of the sensor kit onto the opening of the saline cube.
7. Connect the outlet of the sensor kit with 10mm dia. inlet tube to the pump inlet of the cell washer. Secure the inlet tube with hose band.
8. Connect the sensor cable to the connector at the back of cell washer.

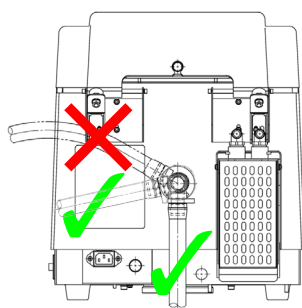


No.	Part
1	Screw cap
2	Inlet tube
3	Liquid sensor
4	Outlet
5	Cord lock
6	Sensor tube
7	Sensor cable

18. Connect waste drain

1. Put the waste drain at a place lower than the cell washer. ②
Make sure that the waste drain is lower than the cell washer to prevent backward flow of waste liquid.
Prepare the saline tank and a drain for waste following your laboratory standards. A waste liquid tank (not included in items supplied) is shown as an example in the picture above.
2. Apply a thin layer of grease (483719) to the inner side of the 18 mm dia. tube (5) or the outer side of the DRAIN connector (4).
3. Connect the 18 mm dia. tube (5) to the DRAIN connector. (5) Use the L-shaped adapter instead if needed.
4. Do not force the 18 mm dia. tube (5) or the L-shaped adapter onto the drain connector (4).
5. Connect the 18 mm dia. tube (5) to the drain connector. ④ Use the L-shaped adaptor instead if needed. Apply a thin layer of grease (483719) to the inner side of the tubing or the outer side of the connector. Do not force the 18 mm dia. tube or the L-shaped adaptor into the drain connector.
Secure the tube or adaptor with the hose band (large).

When the L-shaped adaptor is used, hold it facing downward and connect it to the drain connector. Improper installation may cause waste liquid to stagnate and result in bad drainage.



Make sure to position the L-shaped adapter correctly.

3. 7. Storage

WARNING If necessary clean, disinfect or decontaminate the entire system when removing the cell washer and accessories from use. In doubt contact Thermo Fisher Scientific service.

- Before storing the cell washer and the accessories, it must be cleaned and, if necessary, disinfected and decontaminated.
- The cell washer and accessories have to be thoroughly dried before storage.
- Store the cell washer in a clean and dust-free location.
- Be sure to place the cell washer on its feet.
- Avoid direct sunlight.

3. 8. Shipping

WARNING Before shipping the cell washer and accessories clean and, if necessary, disinfect or decontaminate the entire system. If in doubt, contact Thermo Fisher Scientific service.

Before shipping the cell washer please bear the following in mind:

- The cell washer must be clean and decontaminated.
- The decontamination must be confirmed in a decontamination certificate ([“Decontamination Information Certificate” on page 46](#)). Contact Thermo Fisher Scientific service for more details.
- Install the motor holder for shipping.

4. Operation



WARNING

Use the cell washer only with the rotor, distributor assembly, bowl, splash guard and drain cover installed correctly according to the instructions given in this manual.

Hazardous substances may leak out of the cell washer if the bowl and drain cover are missing.



CAUTION

Open the door carefully after blood cell washing. Remaining waste liquid from the inside of the drain cover may spatter and contaminate persons or the environment.



CAUTION

Close the door by pushing down at the forward center of the door until a short beep is heard.

If the door is not fully closed, an alarm (E1: DOOR OPEN) is indicated and the cell washer does not start. If this alarm is indicated, close the door and press START again.

4. 1. Preparation

1. Prepare the test tubes (10 or 12 mm in diameter and 75 mm in length). When using the 10 mm dia. test tubes, install the D10 adapters (supplied as a standard accessory) to the rotor holder. Make sure that the brim of the adapter is pointing outwards.

CAUTION Be sure to use specified test tubes with a wall thickness of 1mm or more. Test tubes with wall thicknesses of less than 1mm, such as disposable test tubes, are still acceptable to be used as long as the tubes are free from scratches, cracks, warping or other abnormalities that may compromise the structural integrity of the tubes under the centrifugal force exerted by this device, and are not re-used after their initial use is complete. It is recommended to use tempered glass test tube for this cell washer.

2. Check and prepare the saline tank and a drain for waste following your laboratory standards.

CAUTION Make sure that the tips of the 18 mm dia. tubes are above the waste liquid level in the tank. Otherwise, the chamber cannot be drained completely and waste liquid stagnates in the chamber because of poor drainage. Poor drainage can also be caused if the tube is bent or flattened or if waste liquid stagnates in the middle of the tube. Set the tubes properly so that waste liquid drains off smoothly. Be careful not to set the 18 mm dia. tubes at a place higher than the drain outlet of the cell washer.

4. 1. 1. Injection Volume Adjustment and Pump Priming

Prime the Pump

1. With the saline tank filled with buffer saline, hold a beaker or container to the nozzle at the bottom of the lid.



2. Open the cover of the condition-setting panel at the front bottom of the unit.



3. Press SALINE PRIME to discharge saline until the saline coming out of the nozzle is free of bubbles.



NOTICE

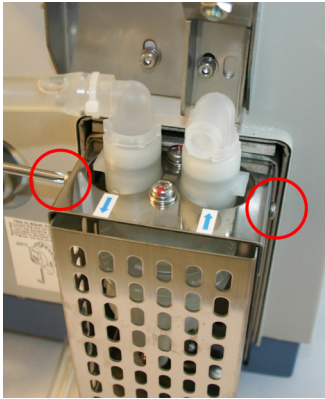
The injection volume is factory set to 96 ml for a 24-place rotor with 12 mm tubes. In case a change of the injection volume is required, please refer to [“4. 1. 1. Injection Volume Adjustment and Pump Priming” on page 34](#).

Set Saline Injection Volume

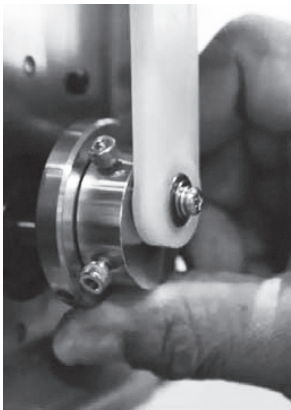
The CW3 cell washer is factory-adjusted to the injection volume of saline for use with the 12 place rotor and the 12 mm diameter test tubes.

When using the 12 place rotor and 12 mm diameter test tubes with this cell washer, there is no need to adjust the injection volume. When using the 24 place rotor or 10 mm diameter test tubes, adjust the flow rate of the pump as follows:

- 1. Turn the power off. Remove the screws as indicated with a screw driver and remove the protective cover of the pump at the back of the device.



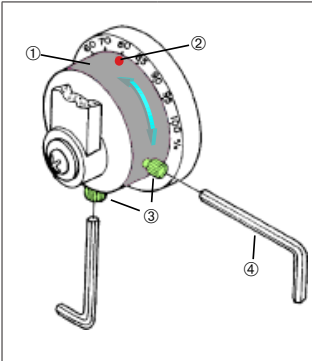
- 2. Loosen the two locking screws on the adjusting knob of the pump anticlockwise, using the hex wrench.



- 3. Turn the adjusting knob to set the indicator value appropriate to your rotor and tube settings, as indicated in the table below.
- 4. Tighten the two locking screws clockwise.

Reference (Index of injection volume)

Rotor	Test tube	Injection volume (+/- 10%)	Indicator setting
12 place	12 mm diameter	48 ml	45 %
12 place	10 mm diameter	32 ml	30 %
24 place	12 mm diameter	96 ml	80 %
24 place	10 mm diameter	64 ml	60 %



The diagram shows a circular pump head assembly. Callout 1 points to the main body of the pump. Callout 2 points to a red dot on a scale around the perimeter. Callout 3 points to two green locking screws on the side. Callout 4 points to a hex wrench used to adjust the screws.

①	Adjusting knob
②	Indicator (red dot)
③	Locking screws
④	Hex wrench

- 5. Saline calibration
 - a. Open the cover of the condition-setting panel at the front bottom of the unit.

- b. Hold a beaker or container against the nozzle at the bottom of the lid and press the OPTION pushbutton. The saline pump runs for 5 seconds.
 - c. Compare the preset volume in the above table and the actual volume injected in the beaker.
 - d. Set the adjusting knob when the actual volume is not in the preset volume range by repeating steps 2 through 4.
 - e. Repeat step a and b until calibration is successful.
6. Reinstall the protective cover of the pump.

Setting the Operating Conditions


The operating conditions are factory-set as follows.

Setting dials and switches	Factory-setting
1 CYCLE	3 cycles
2 AUTO CENTRI TIME	35 seconds
3 MANUAL CENTRI TIME	60 seconds
4 MANUAL SPEED	L (1200 rpm)
5 DECANT SPEED	H-1 (400 rpm)
6 OVERFLOW	1 (Operates only during the first cycle)
7 PROGRAM	1 (5-second centrifugation is performed during the final cycle in automatic mode)
8 AGITATE TIME	5 seconds
9 MELODY	1 (Electronic beep (three beeps))

Agglutination of blood cells is influenced by centrifugal force and centrifugation time. If centrifugal force and centrifugation time are insufficient, agglutination can be insufficient causing incorrect results. Excessive centrifugation can harden the blood cells causing incorrect agglutination or difficult resuspension. The weak reaction may disappear. Centrifugal conditions are influenced by the amount of blood cells, specific gravity of blood cells, etc. The above settings are just a guide. Determine the optimum conditions according to the following evaluation.

1. The supernatant is transparent and no blood cell is suspended after centrifugation.
2. The blood cells that are precipitated at the bottom of the test tubes show clear outlines after centrifugation.
3. The blood cells are easily removed from the bottom and disentangled by a light agitation after decantation.
4. Check the reaction using reagents whether negative or positive.
5. Check the amount of the remaining blood cells. If the amount of the remaining blood cells is small, lower the DECANT SPEED.

4. 1. 2. Preparation of the Sample

	CAUTION
<p>When the sample or liquid spills over, it can leak into the cell washer.</p> <p>If the sample or liquid has leaked into the cell washer, contact Thermo Fisher Scientific service for cleaning and drying.</p>	

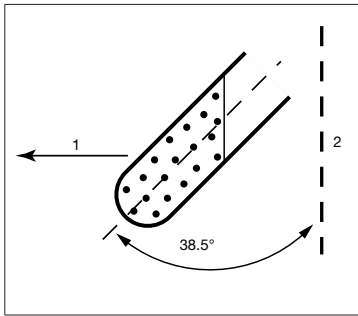
Blood cell washing

Use one or two drops (about 50 µl) of 3-5% erythrocyte suspension in one test tube.

CAUTION When centrifuging precipitated erythrocyte layer, use one or two drops (about 50 µl or less) per test tube. Otherwise, there is a risk that a series of blood cell washing processes cannot be completed.

Centrifugation

Use 80% or less of the test tube capacity as sample volume for centrifugation. The tubes are centrifuged at an angle of 38.5°.



- 1 Centrifugal force
- 2 Drive shaft

4. 2. Operation Modes

4. 2. 1. Automatic Mode

Step	Operation	Description
1	Press the AUTO pushbutton.	AUTO lights up. Automatic mode is selected automatically when the power switch is switched on.
2	Press the START pushbutton.	START lights up. To stop the cell washer temporarily in the middle of a process, press the STOP pushbutton and the cell washer stops at once. For restarting the process, press the START pushbutton.
3	Upon completion of all processes	The end buzzer sounds and START goes out. The door is opened. The buzzer stops as STOP is switched on.

NOTICE In automatic mode, the set number of cycles can be decreased by pushing FEED. For example, if three cycles are set, the number changes 3→2→1 as FEED is held depressed.

4. 2. 2. Manual Mode

In manual mode each process step (WASH, CENTRIFUGE, DECANT, AGITATE) can be run as a single process. More information on the single process steps: „Operational Sequence for Antiglobulin Test“ on page 38.

Step	Operation	Description
1	Switch on MANUAL.	MANUAL lights up.
2	Press FEED once or multiple times to set the desired process manually. The LED of the set PROCESS lights up.	The CENTRIFUGE process is set by default. Possible process steps: <ol style="list-style-type: none"> 1. DECANT - Pressing FEED once (1x) selects the DECANT process. The appropriate LED for the process step lights up. 2. AGITATE - Pressing FEED twice (2x) selects the AGITATE process. The appropriate LED for the process step lights up. 3. WASH - Pressing FEED three times (3x) selects the WASH process. The appropriate LED for the process step lights up. 4. CENTRIFUGE - Pressing FEED four times (4x) selects the CENTRIFUGE process is chosen by default when MANUAL is pressed. The appropriate LED for the process step lights up. Make sure you have set the desired options on the condition-setting panel for the manually controlled process step you are about to launch.
3	Press the START pushbutton.	START lights up.
	Upon completion of the selected process	The end buzzer sounds and the door is opened. START goes out.

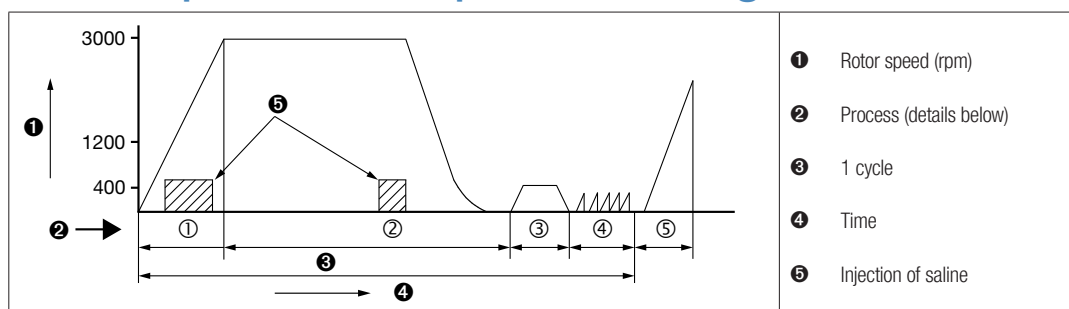
NOTICE The door opens automatically after the rotor is stopped by pressing STOP in the middle of an operation.

NOTICE When power fails during operation and is restored subsequently, or when POWER is switched off during operation and then turned on again, the cell washer stops and reverts to automatic mode operation (AUTO lights up) when power returns. If that happens, set the desired process again and repeat operation in manual mode.

4. 2. 3. Agitation-Centrifugation

Step	Operation	Description
1	Press the AGI/CENT pushbutton.	AGI/CENT lights up.
2	Press the START pushbutton.	START lights up. To stop the cell washer temporarily in the middle of a process, press STOP. For restarting the process, press START.
3	Upon completion of the process	The end buzzer sounds and START goes out. AGI/CENT lights up and the door is opened.

4. 2. 4. Operational Sequence for Antiglobulin Test



	Process	Operation	Description	Reference Illustration
Washing cycle repeated 3 or 4 times	①	WASH	A fixed amount of saline is pumped into the distributor when the rotor speed reaches 1200 rpm. The saline is injected by centrifugal force from the distributor into the test tubes. The blood cells in the test tubes are sufficiently suspended in the saline.	
	②	CENTRIFUGE	Blood cells are centrifuged. The standard centrifugation time is 35 sec. (selectable). Before the rotor decelerates, the injection of saline continues to overflow the test tubes. (Overflow is also selectable.)	
	③	DECANT	The rotor spins at low speed with the rotor holder kept at an angle to slightly open its top end by magnetic force. By this operation only the saline decants from the test tubes and the blood cells remain.	
	④	AGITATE	The rotor repeats rotation and stops at short, quick intervals to disentangle the remaining blood cells.	
	⑤	CENTRIFUGE	The rotor spins about 5 seconds to collect the blood cells adhered to the wall surfaces of the test tubes at the bottom. This is done to ensure the reaction with the Antiglobulin reagent. This operation is performed at the end of the washing cycle.	

5. Maintenance



WARNING

Cell washer, rotor or accessories can be contaminated by samples.
Decontaminate according to good laboratory procedures and methods.



WARNING

If there is a possibility that the cell washer, rotor, or an accessory is contaminated by samples that may impair human health (for example, samples that are toxic or radioactive, or blood samples that are pathogenic or infectious), it is your responsibility to sterilize or decontaminate the cell washer, rotor, or the accessory correctly before requesting repairs from Thermo Fisher Scientific service. Note that Thermo Fisher Scientific can not repair the cell washer, rotor or the accessory unless sterilization or decontamination is completed.



WARNING

It is your responsibility to sterilize and/or decontaminate the cell washer, rotor, or parts correctly before returning them. In such cases copy the decontamination sheet at the end of this manual, fill it out and attach it to the item to be returned. Thermo Fisher Scientific may ask you about the treatment for the cell washer, rotor or the part if the decontamination is checked and judged as insufficient by Thermo Fisher Scientific. It is your responsibility to bear the cost of sterilization or decontamination. Note that Thermo Fisher Scientific can not repair or inspect the cell washer, rotor or the accessory unless sterilization or decontamination is completed.



CAUTION

Do not operate the cell washer in any manner not described in this instruction manual. Should you have any troubles with the cell washer, call Thermo Fisher Scientific service.

5. 1. Cleaning

NOTICE

Information provided is to be considered as general guideline and may vary depending on the usage of the unit.
If you use other cleaning methods than those described here, make sure that the necessary cleanliness is achieved according to your requirements.

5. 1. 1. Pump, Tank and Tubing

CAUTION *If saline is contaminated by bacteria, it may cause hemolysis or poor result.*

- Wash the inside of the tank and the tubing on a regular basis.
- If the cell washer is not used for a long period, drain out saline from the tank, tubing and pump.
- Check the volume of saline inside the tank before operation. Connect the tube securely to the saline tank and the pump inlet connector with the hose bands. When the tube is loose, replace it.
- Make sure that all tubing is properly connected and free from obstructions.

Washing

1. Fill up the 5L saline tank (S413259C) with washing solution (0.5% Sodium Hypochlorite Solution).
2. Fill up the fluid passage with washing solution according to the pump priming procedure outlined in ["Prime the Pump"](#) on page 34.
3. Flush the tubing by running through four wash cycles.



WARNING

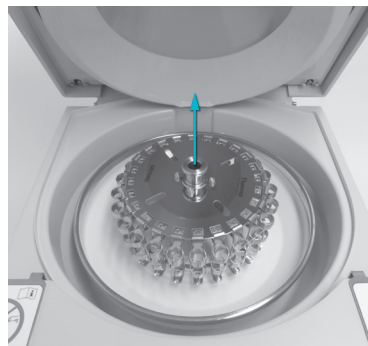
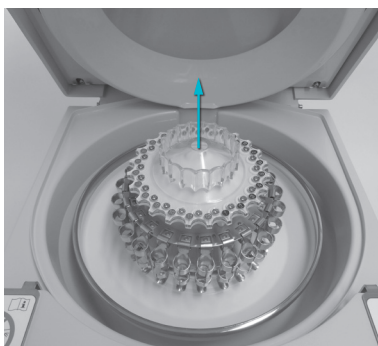
Do not run wash cycles without the rotor and distributor installed. Damage to the bearing may result.

4. Replace the washing solution in saline tank with distilled water
5. Flow 2 to 3 liters of distilled water from the saline tank according to the pump priming procedure outlined in ["Prime the Pump"](#) on page 34.
6. Exchange distilled water in the saline tank with saline.
7. Flow 2 to 3 liters of saline from the saline tank in the same manner as described in step 4.

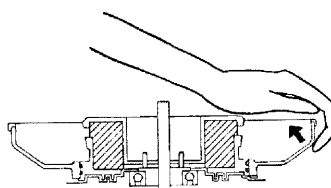
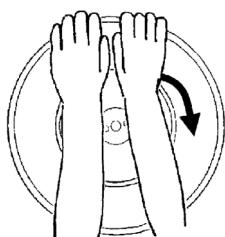
NOTICE *If you are using the Modified saline sensor kit to connect the saline cube with cell washer, disconnect the sensor kit from the cell washer, perform washing steps 1-4. Re-connect the modified saline sensor kit with the cell washer and saline cube. Perform step 6.*

5. 1. 2. Rotor, Distributor and Bowl

After operation remove the rotor, distributor and bowl. Wash and dry them well.



To remove the bowl hold it with both hands. Remove the bowl by lifting it. Slowly rotating the bowl makes it easier to lift.



See the following table for cleaning instructions.

Part	Cleaning
Rotor	Soak in washing solution (0.5% sodium hypochlorite solution). Wash and dry.
Distributor, bowl	Soak in washing solution (0.5% sodium hypochlorite solution) for 1 hour. Wash and dry.

CAUTION The distributor is made of a polycarbonate material. For cleaning make sure that you use washing solution (0.5% sodium hypochlorite solution) to prevent the distributor from deteriorating. Use neutral detergent (pH 6 to 8). Do not soak the distributor in the diluted detergent solution for an extended period of time. Long periods of exposure to detergent may compromise the mechanical strength of the distributor.

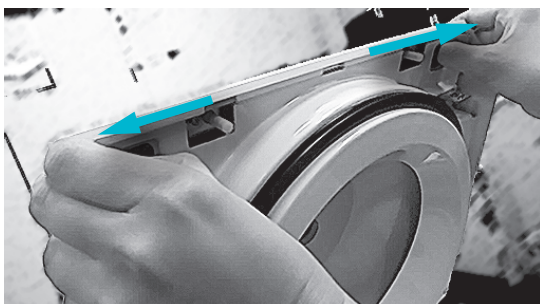
5. 1. 3. Chamber, Splash Guard, Drain Cover and Door Stopper Parts

Remove the splash guard and the drain cover from the cell washer. Wash and dry them well.

Use washing solution (0.5% sodium hypochlorite solution) for cleaning. When washing the inside of the chamber, be careful not to pour washing solution on the decantation coil. Wipe the decantation coil with a cloth or paper moistened with washing solution.

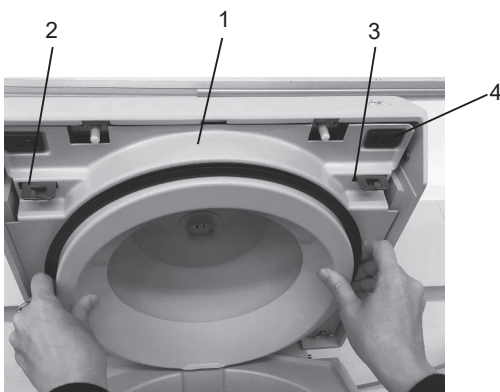
5. 1. 3. 1. Drain Cover

Removing



1. Slide one of the two fasteners to the open position and carefully pull the drain cover out of its seat.
2. Repeat for the second fastener to release the drain cover completely.
3. Pull the drain cover forward to remove it.

Mounting



1. Align the bottom end of the drain cover (1) with the hinges at the rear side of the lid.
2. Align the lid hooks (2) with the slots (3) of the drain cover and push the drain cover firmly into the lid. The drain cover has to be pushed beyond the springs of the hooks.
3. Press the top of the drain cover firmly into place until the two fasteners (4) engage with a clicking sound.

CAUTION When installing the drain cover, make sure that the nozzle at the center of the door goes into the center hole of the drain cover.

5. 1. 3. 2. Splash Guard

Removing

1. Remove the drain cover. See "5. 1. 3. 1. Drain Cover" on page 41.
2. Remove the distributor, rotor and bowl.
3. Remove the splash guard by pulling it up.

Mounting

1. Insert the splash guard.
2. Install the distributor, rotor and bowl. See "3. 6. Set Up" on page 20.
3. Mount the drain cover. See "5. 1. 3. 1. Drain Cover" on page 41.

CAUTION Do not store washing solution in the chamber. Do not spill washing solution on the decantation coil. The liquid may leak into the cell washer causing a failure.

If a test tube is broken, be sure to remove all fragments from inside the chamber and the 18 mm dia. tube of the drain cover thoroughly - for example, using a brush.

5. 1. 3. 3. Checking and Replacing the Door Stopper Parts

If you observe any of the conditions described below, please contact Thermo Fisher Scientific service.

1. The rubber mounts exhibit signs of cracking or other deterioration.
2. The door opens at too wide an angle and remains in that position.

If the door remains wide-open, waste liquid may drip from the drain cover onto the rear panel of the cell washer.

5. 2. Preventive Maintenance

- Replace the 18 mm dia. tube and the 10 mm dia. tube every 1 to 3 years depending on the degree of discoloration.
- The pump is a replacement part. Liquid may start to leak from the bellows of the pump as the pump ages. The degree and speed of deterioration depends on the conditions prevailing in installation environment, such as exposure to ultraviolet rays and high or low temperature. We highly recommended to replace the bellows of the pump (S413230A) every three years.
- Replace the carbon brushes of the motor after 7 years of use (assuming 30 runs per day).

5. 3. Shipping and Disposal



WARNING

When removing the centrifuge and accessories from use for disposal you have to clean and, if necessary, disinfect or decontaminate the entire system. If in doubt contact Thermo Fisher Scientific service.

For the disposal of the centrifuge mind the regulations in your country. Contact Thermo Fisher Scientific service for the disposal of the centrifuge. For contact information check the backpage of this manual or visit www.thermofisher.com/centrifuge

For the countries of the European Union disposal is regulated by the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC.

Observe the information on transport and shipping ("3. Transport and Set Up" on page 19, "3. 8. Shipping" on page 32).

6. Troubleshooting

6. 1. Mechanical Emergency Door Release

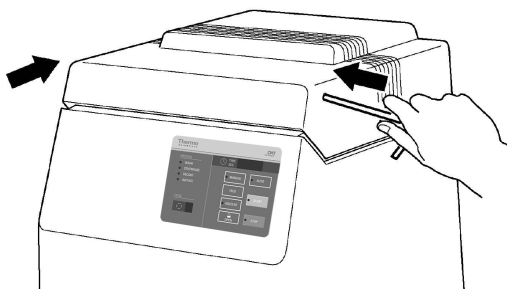


CAUTION

Forcing the door open with the rotor still spinning is an extremely hazardous action. Never unlock the door while the rotor spins. If the door is opened while the rotor still spins, close the door immediately. Do not run the cell washer with the Allen wrench inserted in the hole of the cell washer.

The door cannot be opened and closed except when the cell washer is switched on and the rotor is at a complete standstill. If the door cannot be opened because of a power failure, open the door according to the following procedure.

1. Verify that the rotor has stopped spinning and come to a complete standstill.
Listen carefully to make sure that no spinning sound is heard.
It takes about one minute for a rotor spinning at 3,000 rpm (maximum speed) to stop completely. Make sure you wait long enough before carrying out the next step.
2. To unlock the door, insert the Allen wrench supplied with the unit into the small hole.
A small hole is provided on both sides of the cell washer.
Insert the Allen wrench straight into the small hole and push it in until a clicking sound is heard. Then insert the Allen wrench into the small hole on the other side and repeat the operation. The door unlocks and opens.



6. 2. Error Codes

If any error condition occurs during operation, the corresponding error code blinks on the TIME indicator, the alarm buzzer sounds, and the cell washer stops. Take the action described in the following table or contact Thermo Fisher Scientific service.

Error code	Description	Troubleshooting
E1	The door is open.	Close the door safely and press START.
E2	The saline tank needs to be refilled.	<p>Refill the tank and press STOP to clear the error code. In manual mode, CENTRIFUGE and DECANT processes can be executed even when this error code is shown.</p> <p>If the alarm buzzer sounds even though there is enough saline, the liquid level sensor may be faulty. Contact a Thermo Fisher Scientific authorized sales/service representative for repair. The error code can be temporarily cleared by pressing MANUAL and STOP simultaneously for three seconds or more (operation in automatic mode is possible). It returns to the original state by turning on and off the power switch.</p>
E3	A power failure occurs.	<p>If a power failure occurs while the cell washer is running, the cell washer stops spinning and displays error code "E3".</p> <p>In order to clear the error code, press any of the switches on the control panel after a lapse of one minute or more after the error code has appeared. If the error persists, contact Thermo Fisher Scientific service.</p> <p>This error can be caused by a voltage drop that occurs when the rotor accelerates and your cell washer is connected to a power outlet that serves multiple appliances. Turn off the cell washer power and connect the cell washer to an outlet that can supply stable voltage.</p>
E4	The power supply is incorrect.	The frequency of the power supply is incorrect. Turn off the power switch and on again. If the error persists, contact Thermo Fisher Scientific service.
E5	The rotor overspeeds.	This error code is shown when the rotor runs at a higher than specified speed. Contact Thermo Fisher Scientific service.
E6	An incorrect electric current has been detected.	This error code is shown when an incorrect current flows in the circuitry. Contact Thermo Fisher Scientific service.
E7	The speed sensor is defective.	The sensor for rotor speed is defective. Contact Thermo Fisher Scientific service.
E8	The current sensor is defective.	The sensor for electric current is defective. Contact Thermo Fisher Scientific service.
E9	The RAM is defective.	Turn the power switch off and back on. If the error persists, the microcomputer of the cell washer to control its operation is defective. Contact Thermo Fisher Scientific service.
E10	The rotor overspeeds (hardware detection).	This error code is shown when the rotor runs at a higher than specified speed. Contact Thermo Fisher Scientific service.
E11	The triac is defective.	The electrical component on the circuit board is defective. Contact Thermo Fisher Scientific service.
E14	Set up may be wrong, door lock not engaging.	Check if the rubber gasket and the bowl are installed correctly. „11. Check rubber gasket“ on page 25 and „14. Install bowl“ on page 27. If the error still occurs, contact Thermo Fisher Scientific service.
E16	Set up may be wrong, motor not spinning.	Check if the rubber gasket and the bowl are installed correctly. „11. Check rubber gasket“ on page 25 and „14. Install bowl“ on page 27. If the error still occurs, contact Thermo Fisher Scientific service.
Others	System error	Contact Thermo Fisher Scientific service.

6. 3. Troubleshooting when no Error Code is indicated

	WARNING
Unauthorized disassembly of the cell washer except by Thermo Fisher Scientific service is strictly prohibited.	

No.	Symptom	Possible cause	Remedy
1	POWER does not light its LED even if it is turned on.	1. A power failure occurs. 2. The power cord is disconnected.	1. Wait until power is restored. 2. Contact Thermo Fisher Scientific service.
2	The cell washer does not start operation.	1. The door is not closed. 2. The door switch is defective.	1. Close the door. 2. Contact Thermo Fisher Scientific service.
3	Injection of saline does not take place.	1. No saline is in the tank. 2. The pump is not filled with saline. 3. The pump is defective. 4. The pump connector is loose.	1. / 2. For the first two possible causes: refill the tank with saline to fill the pump. 3. Contact Thermo Fisher Scientific service. 4. Retighten it securely.
4	Saline (waste liquid) scatters in the chamber. 1. The amount of blood cells after washing is less than before. 2. No washing is executed. 3. Washing is executed.	1. Injection volume is excessive and the test tubes are overflowed with solution. 2. The nozzle is out of position and fails to supply the distributor with solution. 3. The drain cover is broken.	1. Adjust the injection volume correctly. 2. / 3. Contact Thermo Fisher Scientific service.
5	No decantation is executed.	The coil is disconnected.	Contact Thermo Fisher Scientific service.
6	Abnormal noise is heard. 1. Contacting noise 2. Others	1. The bowl and the rotor contact the drain cover. 2. The bearing and the locking screw for the decantation coil are defective.	1. Set the bowl and the rotor correctly. 2. Contact Thermo Fisher Scientific service.
7	The cell washer intensely vibrates.	1. Imbalance operation. 2. The bowl, rotor and distributor are not set correctly. 3. The center packing is not mounted correctly. 4. Others.	1. Balance correctly. 2. Set the bowl, rotor and distributor correctly. 3. Mount the center packing correctly. 4. Contact Thermo Fisher Scientific service.
8	A test tube breaks.	1. The test tube is not strong. 2. The test tube is different in size. 3. The rotor holder is deformed.	1. Use test tubes strong enough to bear the centrifugal force. 2. Use applicable-sized test tubes. 3. Contact Thermo Fisher Scientific service.
9	Hemolysis is found in the blood cells.	1. Density of saline is improper. 2. Sample is contaminated by bacteria. 3. Fragments of a test tube are mixed.	1. Exchange for saline of proper density (0.9%). 2. Clean the tank, tube and pump (especially the inside of the tank). 3. Remove glass fragments in the chamber and the drain cover.

Decontamination Information Certificate

INSTRUCTIONS

When an instrument used with radioactive, pathogenic, or otherwise hazardous materials requires servicing by Thermo Fisher Scientific personnel either at the customer's laboratory or at Thermo Fisher Scientific facilities, the following procedure must be complied with to insure safety of our personnel:

1. The instrument or part to be serviced shall be cleaned of all blood and other encrusted material and decontaminated prior to servicing by our representative.
No radioactivity shall be detectable by survey equipment.
2. A Decontamination Information Certificate shall be completed and attached to the instrument or part.

If an instrument or part to be serviced does not have a Decontamination Information Certificate attached to it, and, in our opinion, presents a potential radioactive or biological hazard, our representative will not service the equipment until proper decontamination and certification has been completed.

If an instrument is received at our service facilities and, in our opinion, poses a radioactive or biological hazard, the sender will be contacted for instructions as the equipment is to be disposed. Disposition costs will be borne by the sender.

Additional certificates are available from your local technical or customer service representative. In the event these certificates are not available, a written statement certifying that the instrument or part has been correctly decontaminated and outlining the procedures used will be acceptable.

NOTE *Thermo Fisher Scientific Service representatives will indicate on a Customer Service Repair Report if decontamination was required, and if so, what the contaminate was and what procedure was used. If no decontamination was required, it should be stated.*

Decontamination Information Certificate

Complete and attach to equipment BEFORE servicing.

DECONTAMINATION

CERTIFIED BY _____ TITLE/POSITION _____
 PHONE _____ FAX _____ DEPARTMENT _____
 INSTITUTION ADDRESS _____
 CITY _____ STATE _____ ZIP _____
 INSTRUMENT _____ SERIAL NUMBER _____
 ROTOR _____ SERIAL NUMBER _____
 PART _____ PART NUMBER _____
 HAZARDOUS CONTAMINANT(S) _____ DECONTAMINATION DATE _____
 DECONTAMINATION METHOD(S) _____
 DECONTAMINATION _____
 CERTIFIER'S SIGNATURE: _____ DATE: _____

Chemical Compatibility

CHEMICAL	MATERIAL																											
	ALUMINUM	ANODIC COATING FOR ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	PODIURETHANE ROTOR PAINT	COMPOSITE CARBON FIBER/EPXY	DELTA™	ETHYLENE PROPYLENE	GLASS	NEPRENE	NARL™	NYLON	PET, POLYGLASS™, CLEAR CAMP™	POVALIDMER	POVCARBONATE	POVYESTER, GLASS THERMOSET	POXYTHERMIDE	POXYETHYLENE	POXYPROPYLENE	POXYLUFONE	POXYVINYL CHLORIDE	RLON A™, TELON™	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	Tygon™	Viton™	
2-MERCAPTOETHANOL	S	S	U	/	S	M	S	/	S	U	S	S	U	S	S	/	S	S	S	S	U	S	S	S	S	S	S	
ACETALDEHYDE	S	/	U	U	/	/	/	M	/	U	/	/	/	M	U	U	U	M	M	/	M	S	U	/	S	/	U	
ACETONE	M	S	U	U	S	U	M	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	M	M	S	U	U	
ACETONITRILE	S	S	U	/	S	M	S	/	S	S	U	S	U	M	U	U	/	S	M	U	U	S	S	S	S	U	U	
ALCONOX™	U	U	S	/	S	S	S	/	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S	S	U	
ALYL ALCOHOL	/	/	/	U	/	/	S	/	/	/	/	S	/	S	S	M	S	S	S	/	M	S	/	/	S	/	/	
ALUMINUM CHLORIDE	U	U	S	S	S	S	U	S	S	S	S	M	S	S	S	S	/	S	S	S	S	S	M	U	U	S	S	
FORMIC ACID (100%)	/	S	M	U	/	/	U	/	/	/	/	U	/	S	M	U	U	S	S	/	U	S	/	U	S	/	U	
AMMONIUM ACETATE	S	S	U	/	S	S	S	/	S	S	S	S	S	S	S	U	/	S	S	S	S	S	S	S	S	S	S	
AMMONIUM CARBONATE	M	S	U	S	S	S	S	S	S	S	S	S	S	S	U	U	/	S	S	S	S	S	S	M	S	S	S	
AMMONIUM HYDROXIDE (10%)	U	U	S	U	S	S	M	S	S	S	S	S	/	S	U	M	S	S	S	S	S	S	S	S	S	M	S	
AMMONIUM HYDROXIDE (28%)	U	U	S	U	S	U	M	S	S	S	S	S	U	S	U	M	S	S	S	S	S	S	S	S	S	M	S	
AMMONIUM HYDROXIDE (conc.)	U	U	U	U	S	U	M	S	/	S	/	S	U	S	U	U	S	S	S	/	M	S	S	S	S	/	U	
AMMONIUM PHOSPHATE	U	/	S	/	S	S	S	S	S	S	S	S	/	S	S	M	/	S	S	S	S	S	S	M	S	S	S	
AMMONIUM SULFATE	U	M	S	/	S	S	U	S	S	S	S	S	S	S	S	S	/	S	S	S	S	S	S	U	S	S	U	
AMYL ALCOHOL	S	/	M	U	/	/	S	S	/	M	/	S	/	M	S	S	S	S	M	/	/	/	U	/	S	/	M	
ANILINE	S	S	U	U	S	U	S	M	S	U	U	U	U	U	U	U	/	S	M	U	U	S	S	S	S	U	S	
SODIUM HYDROXIDE (<1%)	U	/	M	S	S	S	/	/	S	M	S	S	/	S	M	M	S	S	S	S	S	S	M	S	S	/	U	
SODIUM HYDROXIDE (10%)	U	/	M	U	/	/	U	/	M	M	S	S	U	S	U	U	S	S	S	S	S	S	M	S	S	/	U	
BARIUM SALTS	M	U	S	/	S	S	S	S	S	S	S	S	S	S	S	M	/	S	S	S	S	S	S	M	S	S	S	
BENZENE	S	S	U	U	S	U	M	U	S	U	U	S	U	U	U	M	U	M	U	U	U	S	U	U	S	U	S	
BENZYL ALCOHOL	S	/	U	U	/	/	M	M	/	M	/	S	U	U	U	U	U	U	/	M	S	M	/	S	/	S	S	
BORIC ACID	U	S	S	M	S	S	U	S	S	S	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	
CESIUM ACETATE	M	/	S	/	S	S	S	/	S	S	S	S	/	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CESIUM BROMIDE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CESIUM CHLORIDE	M	S	S	U	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CESIUM FORMATE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CESIUM IODIDE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CESIUM SULFATE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	
CHLOROFORM	U	U	U	U	S	S	M	U	S	U	U	M	U	M	U	U	U	M	M	U	U	S	U	U	U	M	S	
CHROMIC ACID (10%)	U	/	U	U	S	U	U	/	S	S	S	U	S	S	M	U	M	S	S	U	M	S	M	U	S	S	S	
CHROMIC ACID (50%)	U	/	U	U	/	U	U	/	/	/	/	S	U	U	S	M	U	M	S	S	U	M	S	/	U	M	/	S
CRESOL MIXTURE	S	S	U	/	/	/	S	/	S	U	U	U	U	U	U	/	/	U	U	/	U	S	S	S	S	U	S	
CYCLOHEXANE	S	S	S	/	S	S	S	U	S	U	S	S	U	U	U	M	S	M	U	M	M	S	U	M	M	U	S	
DEIONIZED WATER	S	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	S	S	S	S	
DISTILLED WATER	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	
DEXTRAN	M	S	S	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S	
DIETHYL ETHER	S	S	U	U	S	S	S	U	S	U	U	S	U	U	U	U	U	U	U	U	U	S	S	S	S	M	U	
DIETHYL KETONE	S	/	U	U	/	/	M	/	S	U	/	S	/	M	U	U	U	M	M	/	U	S	/	/	S	U	U	
DIETHYLPYRO-CARBONATE	S	S	U	/	S	S	S	/	S	S	U	S	U	S	U	/	/	S	S	S	M	S	S	S	S	S	S	
DIMETHYLSULFOXIDE	S	S	U	U	S	S	S	/	S	U	S	S	U	S	U	U	/	S	S	U	U	S	S	S	S	U	U	
DIOXANE	M	S	U	U	S	S	M	M	S	U	U	S	U	M	U	U	/	M	M	M	U	S	S	S	S	U	U	
FERRIC CHLORIDE	U	U	S	/	/	/	M	S	/	M	/	S	/	S	/	/	/	S	S	/	/	/	M	U	S	/	S	
ACETIC ACID (GLACIAL)	S	S	U	U	S	S	U	M	S	U	S	U	U	U	U	U	M	S	U	M	U	S	U	U	S	/	U	
ACETIC ACID (5%)	S	S	M	S	S	S	M	S	S	S	S	S	M	S	S	S	S	S	S	S	M	S	S	M	S	S	M	
ACETIC ACID (60%)	S	S	U	U	S	U	S	U	/	S	M	S	U	U	M	U	S	M	S	M	S	M	S	M	U	S	M	U
ETHYL ACETATE	M	M	U	U	S	S	M	M	S	S	U	S	U	M	U	U	/	S	S	U	U	S	M	M	S	U	U	
ETHYL ALCOHOL (50%)	S	S	S	S	S	S	M	S	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	M	S	M	U	
ETHYL ALCOHOL (95%)	S	S	S	U	S	S	M	S	S	S	S	S	U	S	U	/	S	S	S	M	S	S	S	U	S	M	U	
ETHYLENE DICHLORIDE	S	/	U	U	/	/	S	M	/	U	U	S	U	U	U	U	U	U	U	/	U	S	U	/	S	/	S	
ETHYLENE GLYCOL	S	S	S	S	S	S	S	S	S	S	S	S	/	S	U	S	S	S	S	S	S	S	S	M	S	M	S	
ETHYLENE OXIDE VAPOR	S	/	U	/	/	U	/	/	S	U	/	S	/	S	M	/	/	S	S	S	U	S	U	S	S	S	U	
FICOLL-HYPAQUE™	M	S	S	/	S	S	S	/	S	S	S	S	/	S	S	/	S	S	S	S	S	S	S	M	S	S	S	

Troubleshooting when no Error Code is indicated

CHEMICAL	MATERIAL		POLYMER																											
			ALUMINUM	ANODIC COATING FOR ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POURTHANE ROTOR PAINT	COMPOSITE CARBON FIBER/BROV	DELRI™	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NOVEL™	NYLON	PET "PAVECLEAR"™ CLEAR CRAMP™	POLYOLIMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYURETHANE	POLYETHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RUDIN A™ "TELON™"	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON™	VTOR™	
HYDROFLUORIC ACID (10%)	U	U	U	M	/	/	U	/	/	U	U	S	/	S	M	U	S	S	S	S	M	S	U	U	U	/	/			
HYDROFLUORIC ACID (50%)	U	U	U	U	/	/	U	/	/	U	U	U	U	S	U	U	U	S	S	M	M	S	U	U	U	/	M			
HYDROCHLORIC ACID (30%)	U	U	U	U	/	U	U	M	/	U	M	U	U	M	U	U	U	/	S	/	U	S	U	U	U	/	/			
FORMALDEHYDE (40%)	M	M	M	S	S	S	S	M	S	S	S	S	M	S	S	S	U	S	S	M	S	S	S	M	S	M	U			
GLUTARALDEHYDE	S	S	S	S	/	/	S	/	S	S	S	S	S	S	S	/	/	S	S	S	/	/	S	S	S	/	/			
GLYCEROL	M	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S			
GUMMINE HYDROCHLORIDE	U	U	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	U	S	S	S			
HAEMO-SOL™	S	S	S	/	/	/	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	S	S	S	S			
HEXANE	S	S	S	/	S	S	S	/	S	S	U	S	U	M	U	S	S	U	S	S	M	S	U	S	S	U	S			
ISOBUTYL ALCOHOL	/	/	M	U	/	/	S	S	/	U	/	S	U	S	S	M	S	S	S	/	S	S	S	/	S	/	S			
ISOPROPYL ALCOHOL	M	M	M	U	S	S	S	S	S	U	S	S	U	S	U	M	S	S	S	S	S	S	S	M	M	M	S			
IODOACETIC ACID	S	S	M	/	S	S	S	/	S	M	S	S	M	S	S	/	M	S	S	S	S	S	M	S	S	M	M			
POTASSIUM BROMIDE	U	S	S	/	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S	/	S	S	S	M	S	S	S			
POTASSIUM CARBONATE	M	U	S	S	S	S	S	/	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S			
POTASSIUM CHLORIDE	U	S	S	/	S	S	S	S	S	S	S	S	S	S	S	/	S	S	S	S	S	S	S	U	S	S	S			
POTASSIUM HYDROXIDE (5%)	U	U	S	S	S	S	M	/	S	S	S	S	/	S	U	S	S	S	S	S	S	S	M	U	M	S	U			
POTASSIUM HYDROXIDE (30%)	U	U	M	U	/	/	M	/	M	S	S	/	U	M	U	U	U	S	M	/	M	U	/	U	U	/	U			
POTASSIUM PERMANGANATE	S	S	S	/	S	S	S	/	S	S	S	U	S	S	S	M	/	S	M	S	U	S	S	M	S	U	S			
CALCIUM CHLORIDE	M	U	S	S	S	S	S	S	S	S	S	S	S	S	M	S	/	S	S	S	S	S	M	S	S	S	S			
CALCIUM HYPOCHLORITE	M	/	U	/	S	M	M	S	/	M	/	S	/	S	M	S	/	S	S	S	M	S	M	U	S	/	S			
KEROSENE	S	S	S	/	S	S	S	U	S	M	U	S	U	M	M	S	/	M	M	M	S	S	U	S	U	S	S			
SODIUM CHLORIDE (10%)	S	/	S	S	S	S	S	S	/	/	/	S	S	S	S	S	/	S	S	S	S	/	S	S	M	/	S			
SODIUM CHLORIDE (54%o)	U	/	S	U	S	S	S	/	/	/	/	S	S	S	S	S	/	S	S	/	S	/	S	S	M	/	S			
CARBON TETRACHLORIDE	U	U	M	S	S	U	M	U	S	U	U	S	U	M	U	S	S	M	M	S	M	M	M	M	U	S	S			
AQUA REGIA	U	/	U	U	/	/	U	/	/	/	/	/	U	U	U	U	U	U	U	/	/	/	/	/	S	/	M			
SOLUTION 555 (20%)	S	S	S	/	/	/	S	/	S	S	S	S	S	S	S	/	/	S	S	S	/	S	S	S	S	S	S			
MAGNESIUM CHLORIDE	M	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S	S			
MERCAPTOACETIC ACID	U	S	U	/	S	M	S	/	S	M	S	U	U	U	U	/	S	U	U	S	M	S	U	S	S	S	S			
METHYL ALCOHOL	S	S	S	U	S	S	M	S	S	S	S	S	U	S	U	M	S	S	S	S	S	S	M	S	M	U	U			
METHYLENE CHLORIDE	U	U	U	U	M	S	S	U	S	U	U	S	U	U	U	U	U	M	U	U	U	S	S	M	U	S	U			
METHYL ETHYL KETONE	S	S	U	U	S	S	M	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	S	S	S	U	U			
METRAMIDE™	M	S	S	/	S	S	S	/	S	S	S	S	/	S	S	/	/	S	S	S	S	S	S	M	S	S	S			
LACTIC ACID (100%)	/	/	S	/	/	/	/	/	/	M	S	U	/	S	S	S	M	S	S	/	M	S	M	S	S	/	S			
LACTIC ACID (20%)	/	/	S	S	/	/	/	/	/	M	S	M	/	S	S	S	S	S	S	M	S	M	S	S	S	/	S			
N-BUTYL ALCOHOL	S	/	S	U	/	/	S	/	/	S	M	/	U	S	M	S	S	S	S	M	M	S	M	/	S	/	S			
N-BUTYL PHTHALATE	S	S	U	/	S	S	S	/	S	U	U	S	U	U	U	M	/	U	U	S	U	S	M	M	S	U	S			
N, N-DIMETHYLFORMAMIDE	S	S	S	U	S	M	S	/	S	S	U	S	U	S	U	U	/	S	S	U	U	S	M	S	S	S	U			
SODIUM BORATE	M	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S			
SODIUM BROMIDE	U	S	S	/	S	S	S	/	S	S	S	S	S	S	S	S	/	S	S	S	S	S	S	M	S	S	S			
SODIUM CARBONATE (2%)	M	U	S	S	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S			
SODIUM DODECYL SULFATE	S	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S			
SODIUM HYPOCHLORITE (5%)	U	U	M	S	S	M	U	S	S	M	S	S	S	M	S	S	S	S	M	S	S	S	M	U	S	M	S			
SODIUM IODIDE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S			
SODIUM NITRATE	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	/	S	S	S	S	S	U	S	S	S	S			
SODIUM SULFATE	U	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S	S			
SODIUM SULFIDE	S	/	S	S	/	/	/	S	/	/	/	S	S	S	U	U	/	/	S	/	/	/	S	S	M	/	S			
SODIUM SULFITE	S	S	S	/	S	S	S	S	M	S	S	S	S	S	S	M	/	S	S	S	S	S	S	S	S	S	S			
NICKEL SHOTS	U	S	S	S	S	S	/	S	S	S	/	/	S	S	S	S	/	S	S	S	S	S	S	M	S	S	S			
OILS (PETROLEUM)	S	S	S	/	/	/	S	U	S	S	S	S	U	U	M	S	M	U	U	S	S	S	U	S	S	S	S			
OILS (OTHER)	S	/	S	/	/	/	S	M	S	S	S	S	U	S	S	S	S	U	S	S	S	S	/	S	S	M	S			
OLEIC ACID	S	/	U	S	S	S	U	U	S	U	S	S	M	S	S	S	S	S	S	S	S	S	M	U	S	M	M			
OXALIC ACID	U	U	M	S	S	S	U	S	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	U	M	S	S			
PERCHLORIC ACID (10%)	U	/	U	/	S	U	U	/	S	M	M	/	/	M	U	M	S	M	M	/	M	S	U	/	S	/	S			
PERCHLORIC ACID (70%)	U	U	U	/	/	U	U	/	S	U	M	U	U	M	U	U	U	M	M	U	M	S	U	U	S	U	S			
PHENOL (5%)	U	S	U	/	S	M	M	/	S	U	M	U	U	S	U	M	S	M	S	U	U	S	U	M	M	M	S			
PHENOL (50%)	U	S	U	/	S	U	M	/	S	U	M	U	U	U	U	U	S	U	M	U	U	S	U	U	U	M	S			

CHEMICAL	MATERIAL																													
	ALUMINUM	ANODIC COATING FOR ALUMINUM		BUNA N	CELLULOSE ACETATE BUTIRATE		POLYURETHANE ROTCH PAINT		COMPOSITE CARBON FIBER/EPoxy		DELTA™	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NOPL™	NYLON	PET*, POLYGLASS™, CLEAR CRAMP™	POVALOMER	POLYCARBONATE	POLYESTER GLASS THERMOSET	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	POLYURETHANE	
PHOSPHORIC ACID (10%)	U	U	M	S	S	S	S	U	S	S	S	S	U	/	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S
PHOSPHORIC ACID (CONC.)	U	U	M	M	/	/	U	S	/	M	S	U	U	M	M	S	S	S	S	M	S	M	S	U	M	U	/	S		
PARALOG: MED (SERUM UREA)	M	S	S	S	/	/	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
PICRIC ACID	S	S	U	/	S	M	S	S	S	M	S	U	S	S	S	U	S	S	S	S	S	S	S	U	S	U	M	S	M	S
PYRIDINE (50%)	U	S	U	U	S	U	U	/	U	S	S	U	U	M	U	U	/	U	S	M	U	S	S	U	U	U	U	U	U	U
RUBIDIUM BROMIDE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	S	M	S	S	S	S
RUBIDIUM CHLORIDE	M	S	S	/	S	S	S	/	S	S	S	S	S	S	S	S	/	/	S	S	S	S	S	S	M	S	S	S	S	S
SUCROSE	M	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
SUCROSE, ALKALINE	M	S	S	/	S	S	S	/	S	S	S	U	S	S	U	S	S	S	S	S	S	S	S	S	M	S	S	S	S	S
SULFOSALICYLIC ACID	U	U	S	S	S	S	S	/	S	S	S	U	S	S	S	/	S	S	/	S	S	/	S	S	U	S	U	S	S	S
NITRIC ACID (10%)	U	S	U	S	S	U	U	/	S	U	S	U	/	S	S	S	S	S	S	S	S	S	S	M	S	S	S	S	S	S
NITRIC ACID (50%)	U	S	U	M	S	U	U	/	S	U	S	U	U	M	M	U	M	M	M	S	S	S	S	U	S	S	M	S	S	S
NITRIC ACID (95%)	U	/	U	U	/	U	U	/	/	U	U	U	U	M	U	U	U	U	M	U	U	M	U	S	S	/	S			
HYDROCHLORIC ACID (10%)	U	U	M	S	S	S	U	/	S	S	S	U	U	S	U	S	S	S	S	S	S	S	S	S	U	M	S	S	S	S
HYDROCHLORIC ACID (50%)	U	U	U	U	S	U	U	/	S	M	S	U	U	M	U	U	S	S	S	S	S	S	M	S	M	U	U	M	M	S
SULFURIC ACID (10%)	M	U	U	S	S	U	U	/	S	S	M	U	S	S	S	S	S	S	S	S	S	S	S	S	U	U	U	S	S	S
SULFURIC ACID (50%)	M	U	U	U	S	U	U	/	S	S	M	U	U	S	U	U	M	S	S	S	S	S	S	U	U	U	M	S	S	S
SULFURIC ACID (CONC.)	M	U	U	U	/	U	U	M	/	/	M	U	U	S	U	U	U	M	S	U	M	S	U	U	U	/	S			
STEARIC ACID	S	/	S	/	/	/	S	M	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S	M	M	S	S	S	S	S
TETRAHYDROFURAN	S	S	U	U	S	U	U	M	S	U	U	S	U	U	U	/	M	U	U	U	U	U	S	U	S	S	U	U	U	U
TOLUENE	S	S	U	U	S	S	M	U	S	U	U	S	U	U	U	S	U	M	U	U	U	U	S	U	S	U	U	M	S	S
TRICHLOROACETIC ACID	U	U	U	/	S	S	U	M	S	U	S	U	U	S	M	/	M	S	S	U	U	S	U	U	U	M	U	U	U	U
TRICHLOROETHANE	S	/	U	/	/	/	M	U	/	U	/	S	U	U	U	U	U	U	U	U	U	U	S	U	/	S	/	S	S	
TRICHLOROETHYLENE	/	/	U	U	/	/	/	U	/	U	/	S	U	U	U	U	U	U	U	U	U	U	S	U	/	U	/	S	S	
TRISODIUM PHOSPHATE	/	/	/	S	/	/	M	/	/	/	/	/	/	S	/	/	S	S	S	/	/	S	/	/	S	/	S	/	S	S
TRIS BUFFER (NEUTRAL pH)	U	S	S	S	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
TRITON X/100™	S	S	S	/	S	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
UREA	S	/	U	S	S	S	S	/	/	/	/	S	S	S	M	S	S	S	S	/	S	S	S	M	S	/	S			
HYDROGEN PEROXIDE (10%)	U	U	M	S	S	U	U	/	S	S	S	U	S	S	M	U	S	S	S	S	S	S	S	S	M	S	U	S	S	S
HYDROGEN PEROXIDE (3%)	S	M	S	S	S	/	S	/	S	S	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S	S	S	S
XYLENE	S	S	U	S	S	S	M	U	S	U	U	U	U	U	U	M	U	M	U	U	U	U	S	U	M	S	U	S	S	S
ZINC CHLORIDE	U	U	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	S
ZINC SULFATE	U	S	S	/	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
CITRIC ACID (10%)	M	S	S	M	S	S	M	S	S	S	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S	S	S	S

* Polyethyleneterephthalate

S – Satisfactory.

M – Moderate attack, may be satisfactory for use in centrifuge depending on length of exposure, speed involved, etc.; suggest testing under actual conditions of use.

U – Unsatisfactory, not recommended.

/ – Performance unknown; suggest testing, using sample to avoid loss of valuable material.

NOTICE

Chemical resistance data is included only as a guide to product use. Because no organized chemical compatibility data exists for materials under the stress of centrifugation, when in doubt we recommend pretesting sample lots.

Index

A

Agitation-Centrifugation 38
Antiglobulin Test 38
Automatic Mode 37

B

Before Setting Up 19
Bowl 40

C

Chamber 41
Chemical Compatibility 47
Cleaning 39
Condition-setting Panel 17
Control Panel 16

D

Decontamination Information Certificate 46
Directives and Standards 13
Disposal 42
Distributor 40
Door Stopper 42
Door Stopper Parts 41
Drain Cover 41

E

Error Codes 44

I

Injection Volume Adjustment 34
Intended Use 7
Items Supplied 5

L

Leveling 20
Location 19
Location and Function of Parts 14

M

Mains Connection 20
Mains Supply 13
Maintenance 39
Manual Mode 37
Mechanical Emergency Door Release 43

O

Operating Conditions 36
Operation 33
Operation Modes 37

P

Precautions 7
Preface 5
Preparation 33
Preparation of the Sample 36
Pump 40
Pump Priming 34

R

Rotor 40

S

Setting Up 20
Set Up 19
Shipping 32, 42
Splash Guard 41, 42
Storage 32
Symbols Used in the Manual 11
Symbols Used on the Cell Washer 10

T

Tank 40
Technical Specifications 12
Transport 19
Transporting 19
Troubleshooting 43
Troubleshooting 45
Tubing 40



Thermo Electron LED GmbH
Zweigniederlassung Osterode
Am Kalkberg, 37520 Osterode am Harz
Germany



thermofisher.com/centrifuge

© 2015-2021 Thermo Fisher Scientific Inc. All rights reserved.

Delrin, TEFLON, and Viton are registered trademarks of DuPont. Noryl is a registered trademark of SABIC. POLYCLEAR is a registered trademark of Hongye CO., Ltd. Hypaque is a registered trademark of Amersham Health As. RULON A and Tygon are registered trademarks of Saint-Gobain Performance Plastics. Alconox is a registered trademark of Alconox. Ficoll is a registered trademark of GE Healthcare. Haemo-Sol is a registered trademark of Haemo-Sol. Triton X-100 is a registered trademark of Sigma-Aldrich Co. LLC. Valox is a registered trademark of General Electric Co.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details. Shown pictures within the manual are examples and may differ considering the set parameters and language.

United States/Canada +1 866 984 3766
Latin America +1 866 984 3766
Austria +43 1 801 40 0
Belgium +32 53 73 42 41
France +33 2 2803 2180
Germany national toll free 0800 1 536 376
Germany international +49 61 84 90 6000
Italy +39 02 95059 552

Netherlands +31 76 579 55 55
Nordic/Baltic States/CIS +358 10 329 2200
Russia +7 812 703 42 15
Spain/Portugal +34 93 223 09 18
Switzerland +41 44 454 12 22
UK/Ireland +44 870 609 9203
India +91 22 6716 2200
China +800 810 5118 or

+400 650 5118
Japan +81 3 5826 1616
Other Asian Countries +852 2885 4613
Australia +61 39757 4300
New Zealand +64 9 980 6700
Countries not listed +49 6184 90 6000

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