Thermo Scientific Sorvall Legend X1R Centrifuge

Instruction Manual

50156051_b • 03 / 2020



Visit us online to register your warranty: thermofisher.com/labwarranty

Table of Contents

	Preface	iii
	Scope of supply	iii
	Intended use	iv
	Accident prevention	iv
	Precautions	v
Chapter 1	Introduction and description	1-1
-	Characteristics of the Thermo Scientific Sorvall Legend X1R	
	Technical data	
	Directives, standards and guidelines	
	Functions and features	
	Mains supply	
	Rotor selection	1-6
Chapter 2	Before use	
	Before setting up	
	Transporting the centrifuge	
	Location	
	Aligning the centrifuge	
	Mains Connection	
	Storage	
Chapter 3	Control panel	
	Control panel	
	Keys	
Chapter 4	Operation	4-1
•	Switch on centrifuge	
	Door opening	
	Close door	
	Rotor installation	
	Entering parameters	
	Centrifugation	
	Temperature adaptation during standstill	
	Short-term centrifugation	
	Removing the rotor	
	Aerosol-tight rotors	
	Turning off the centrifuge	
	00-	

1

Chapter 5	System menu	5-1
•	System menu Description	
Chapter 6	Maintenance and care	6-1
-		
	Cleaning intervals	
	Cleaning	
	Disinfection	
	Decontamination	
	Autoclaving	
	Service of Thermo Fisher Scientific	
Chapter 7	Troubleshooting	
	Mechanical emergency lid release	
	Troubleshooting by user	
	When to contact customer service	
	Chemical compatibility chart	A-1
	Contact information	B-1

Preface

Before starting to use the centrifuge, read through this instruction manual carefully and follow the instructions.

The information contained in this instruction manual is the property of Thermo Fisher Scientific; it is forbidden to copy or pass on this information without explicit approval.

Failure to follow the instructions and safety information in this instruction manual will result in the expiration of the sellers warranty.

Scope of supply

Article number		Quantity	Check
	Sorvall Legend X1R Centrifuge	1	
	Power supply cable	1	
50120979	Instruction manual	1	
50122054	CD with manual	1	

If any parts are missing, please contact your nearest Thermo Fisher Scientific representative.



This symbol refers to general hazards.

CAUTION means that material damage could occur.

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



This symbol refers to biological hazards.

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

Intended use

• This centrifuge is designed to separate sample mixtures of different densities like chemicals, environmental samples and other non-human body samples.

Accident prevention



WARNING Plug the centrifuge only into sockets which have been properly grounded.



WARNING If a hazardous situation occurs, turn off the power supply to the centrifuge and leave the area immediately.

Prerequisite for the safe operation of the Sorvall Legend X1R is a work environment in compliance with standards, directives and trade association safety regulations and proper instruction of the user.

The safety regulations contain the following basic recommendations:



- Maintain a radius of at least 30 cm around the centrifuge.
- Implementation of special measures which ensure that no one can approach the centrifuge for longer than absolutely necessary while it is running.

The mains plug must be freely accessible at all times. Pull out the power supply plug or disconnect the power supply in an emergency.

Precautions

In order to ensure safe operation of the Sorvall Legend X1R, the following general safety regulations must be followed:

- The centrifuge should be operated by trained specialists only.
- The centrifuge is to be used for its intended use only.
- Do not move the centrifuge while it is running.
- Do not lean on the centrifuge.
- Use only rotors and accessories for this centrifuge which have been approved by Thermo Fisher Scientific. Exceptions to this rule are commercially available glass or plastic centrifuge tubes, provided they have been approved for the speed or the RCF value of the rotor.
- Do not use rotors which show any signs of corrosion and/or cracks.
- Do not touch the mechanical components of the rotor and do not make any changes to the mechanical components.
- Use only with rotors which have been properly installed. Follow the instructions on the Auto-Lock in section "Rotor installation" on page 4-2.



- Use only with rotors which have been loaded properly. Follow the instructions given in the rotor manual.
- Never overload the rotor. Follow the instructions given in the rotor manual.
- Never start the centrifuge when the lid is open.
- Never open the lid until the rotor has come to a complete stop and this has been confirmed in the display.
- The lid emergency release may be used in emergencies only to recover the samples from the centrifuge, e.g. during a power failure (see section "Mechanical emergency lid release" on page 7-2).
- Never use the centrifuge if parts of its cover panels are damaged or missing.
- Do not touch the electronic components of the centrifuge or alter any electronic or mechanical components.
- Please observe the safety instructions.

Please pay particular attention to the following aspects:

- Location: well-ventilated environment, set-up on a level and rigid surface with adequate load-bearing capacity.
- Rotor installation: Make sure the rotor is locked properly into place before operating the centrifuge.

v



- Especially when working with corrosive samples (salt solutions, acids, bases), the accessory parts and vessel have to be cleaned carefully.
- Always balance the samples.

Centrifuging hazardous substances:

- Do not centrifuge explosive or flammable materials or substances which could react violently with one another.
- The centrifuge is neither inert nor protected against explosion. Never use the centrifuge in an explosion-prone environment.
- Do not centrifuge inflammable substances.

Remaining risk: Improper use can cause damages, contamination, and injuries with fatal consequences.



• Do not centrifuge toxic or radioactive materials or any pathogenic micro-organisms without suitable safety precautions.

When centrifuging microbiological samples from the Risk Group II (according to the Bio-safety Manual" of the World Health Organization (WHO)), aerosol-tight biological seals have to be used.

For materials in a higher risk group, extra safety measures have to be taken.

• If toxins or pathogenic substances have gotten into the centrifuge or its parts, appropriate disinfection measures have to be taken (see "Disinfection" on page 6-4).

Remaining risk: Improper use can cause damages, contamination, and injuries with fatal consequences.

• Highly corrosive substances which can cause material damage and impair the mechanical stability of the rotor, should only be centrifuged in corresponding protective tubes.

Introduction and description

Contents

- "Characteristics of the Thermo Scientific Sorvall Legend X1/ X 1R" on page 1-2
- "Technical data" on page 1-3
- "Directives, standards and guidelines" on page 1-4
- "Functions and features" on page 1-5
- "Mains supply" on page 1-5
- "Rotor selection" on page 1-6

1

Characteristics of the Thermo Scientific Sorvall Legend X1R

Several rotors with a wide range of tubes can be used.

The set speed is reached within seconds. The maintenance-free induction motor ensures quiet and low-vibration operation even at high speeds, and guarantees a very long lifetime.

The user-friendly control panel makes it easy to pre-set the speed, RCF value, run time, temperature, and run profile (acceleration and braking curves). You can choose between the display of speed and RCF or the entry mode.

These settings can be changed even while the centrifuge is running.

With the help of the 🚵 key , you can also centrifuge a sample for just a few seconds, if called for.

The Sorvall Legend X1R is equipped with various safety features:

- The housing and rotor chamber consist of steel plate, the interior of armour steel, while the front panel is made of high-impact resistant plastic.
- The lid is equipped with a view port and a lock.
- The lid of the centrifuge can only be opened while the centrifuge is switched on and the rotor has come to a complete stop. The centrifuge cannot be started until the lid has been closed properly.
- The integrated rotor detection systems ensures that no inadmissible speed settings can be preselected.
- Electronic imbalance recognition
- Lid emergency release: For emergencies only, e.g. during power failures (see "Mechanical emergency lid release" on page 7-2)

Technical data

The technical data of the Sorvall Legend X1R is listed in the following table. **Table 1-1.** Technical Data Sorvall Legend X1R

Feature	Sorvall Legend X1R
Environmental conditions	-Use in interior spaces
	-Altitudes of up to 2,000m above sea level
	-max. relative humidity 80% up to 31 °C; decreasing linearly up to 50% relative humidity at 40 °C.
Permissible ambient temperature	+2 °C to +35 °C
Overvoltage category	II
Pollution degree	2
Voltage	230 V
Heat dissipation	4778 BTU/h
10	
IP	20
Run time	unlimited
Max speed n _{max}	15200 rpm (depending on the rotor)
Min speed n _{min}	300 rpm
Maximum RCF value at n _{max}	25830 x g
Maximum kinetic energy	62.5 kJ
Noise level at maximum speed	< 63 dB (A)
Temperature setting range	-10 °C to +40 °C
Dimensions	refrigerated
Height	360 mm
Height with lid open	870 mm
Width	623 mm
Depth	605 mm
Table top height	310 mm
Weight without rotor	91.5 kg

Refrigerant

Order No.	Centrifuge	Refrigerant	Quantity	Pressure	GWP CO2e
75004260	Thermo Scientific Sorvall Legend X1R	R-404A	0.38 kg	31 bar	3922 1.49 t

Contains fluorinated greenhouse gases in a hermetically sealed system.

Directives, standards and guidelines

Table 1-2. Directives, standards and guidelines

Tension / frequency	Standards
220 / 230 V 50 / 60 Hz	IEC 61010-1 IEC 61010-2-020
	EN 61326-1 Class B
	EN ISO 9001

Functions and features

The following table gives an overview of the important functional and performance characteristics of the Sorvall Legend X1R.

Table 1-3. Functions and Features

Component / function	Description / features
Structure / Housing	Galvanized steel chassis with armoured plating
Chamber	Stainless steel
Drive	Induction drive without carbon brushes
Keys and display	Easy-to-clean keypad and display surface
Controls	Microprocessor-controlled
Internal memory	The most recent data is saved
Functions	RCF-selection, temperature control, pretemp with cooling equipped devices
Acceleration / braking profiles	9 acceleration and 10 braking curves
Rotor recognition	Automatic
Imbalance recognition	Electronic, contingent on rotor and speed
Lid lock	Automatic lid closing and locking starting from an initial hold position

Mains supply

The following table contains an overview of the electrical connection data for the Sorvall Legend X1R. This data is to be taken into consideration when selecting the mains connection socket.

Table 1-4. Electrical connection data of the Sorvall Legend X1R

Cat.	Mains voltage	Frequency	Rated current	Power consumption	Equipment fuse	Building fuse
refrigerated	230 V ±10 %	50 / 60 Hz	8 A	1400 W	15 AT	16 AT

¹Please contact the Thermo Fisher service to run the unit at 60Hz Line frequency.

Rotor selection

The Sorvall Legend X1R of the is supplied without a rotor.

Various Thermo Scientific rotors are available to choose from.

TX-400	7500 3629
with round buckets	7500 3655
TX-200	7500 3658
with round buckets	7500 3659
BIOShield TM 720	7500 3621
M-20	7500 3624
Microliter 30x2 sealed	7500 3652
CLINIConic [™] 30x15	7500 3623
8x50 sealed	7500 3694
Fiberlite [®] F15-6x100	7500 3698
Fiberlite F13-14x50c	7500 3661
Fiberlite F15-8x50c	7500 3663
Fiberlite F21-48x2	7500 3664
Microliter 48x2 sealed	7500 3602
HIGHConic TM II	7500 3620

The technical data of the rotors and the corresponding adapters and reduction sleeves for various commercially available containers can be found in the corresponding rotor operating manuals.

For more information visit our website at: http://www.thermofisher.com/centrifuge

Before use

Contents

- "Before setting up" on page 2-2
- "Transporting the centrifuge" on page 2-2
- "Location" on page 2-2
- "Aligning the centrifuge" on page 2-3
- "Mains Connection" on page 2-3
- "Storage" on page 2-4

2

Before setting up

- Check the centrifuge 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 the order for completeness (see "Scope of supply" on page iii). If the order is incomplete, please contact Thermo Fisher Scientific.

Transporting the centrifuge

- Due to its weight (see "Technical data" on page 1-3), the centrifuge should be carried by several people.
- Always lift the centrifuge at both sides.



Figure 2-1. Lifting the centrifuge at both sides.

- The centrifuge can be damaged by impacts.
- Transport the centrifuge upright and if at all possible in its packaging.



WARNING Always lift the centrifuge on both sides. Never lift the centrifuge by its front or the back panel. Always remove the rotor before moving the centrifuge.

Location

The centrifuge should only be operated indoors.

The set-up location must fullfil the following requirements:

- A safety zone of at least 30 cm must be maintained around the centrifuge. People and hazardous substances must be kept out of the safety zone while centrifuging.
- The supporting structure must be stable and free of resonance, for example a level laboratory bench.
- The supporting structure must be suitable for horizontal setup of the centrifuge.

• The centrifuge should not be exposed to heat and strong sunlight.



WARNING UV rays reduce the stability of plastics. Do not subject the centrifuge, rotors and plastic accessories to direct sunlight.

• The set-up location must be well-ventilated at all times.

Aligning the centrifuge

The horizontal alignment of the centrifuge must be checked every time after moving it to a different location.

The supporting structure must be suitable for horizontal setup of the centrifuge.



CAUTION If the centrifuge isn't level, imbalances can occur and the centrifuge can be damaged.

Do not place anything under the feet to level the centrifuge.

Mains Connection

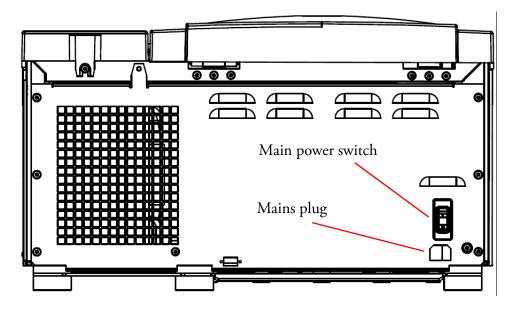


Figure 2-2. Mains connection

- 1. Turn off the power supply switch on the back (press "0").
- 2. Plug the centrifuge into grounded electrical sockets only.
- 3. Check whether the cable complies with the safety standards of your country.
- 4. Make sure that the voltage and frequency correspond to the figures on the rating plate.
- 5. Establish the connection to the power supply with the connecting cable.

Storage

- Before storing the centrifuge and the accessories it must be cleaned and if necessary disinfected and decontaminated.
- Store the centrifuge in a clean, dust-free location.
- Be sure to place the centrifuge on its feet.
- Avoid direct sunlight.

Control panel

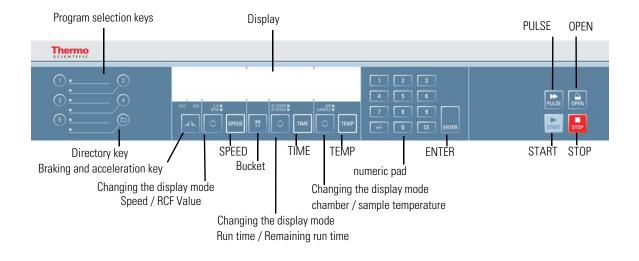
Contents

- "Control panel" on page 3-2
- "Keys" on page 3-2

3

Control panel

The control panel contains the keys and displays of the centrifuge (only the power switch is located on the back of the device). All parameters can be selected and changed during operation.



Keys

The keys allow user intervention for controlling the operating mode as follows:

Key		Display contents
START	Start	Normal start of the centrifuge
STOP	Stop	End run manually
OPEN	Open lid	Automatic release (possible only when device is switched on); Emergency release (see "Mechanical emergency lid release" on page 7-2)
PULSE	Pulse	By pressing the 🚵 key the centrifuge starts immediately and accelerates up to the end speed. Releasing the key initiates a stopping process at the highest braking curves.
	Directory	By pressing the 🕞 key up and down you change the value in thedisplay.
T	Bucket	Use the key in order to have all available bucket types displayed in succession.
	Braking and accelera- tion curves	Use the 🔼 key to toggle between braking and acceleration curves.
	Changing the display mode	Use the key 这 to change the display mode. (speed / RCF value, sample / chamber temperature, run time counter from start or preset speed on)

Operation

Contents

- "Switch on centrifuge" on page 4-2
- "Door opening" on page 4-2
- "Rotor installation" on page 4-2
- "Entering parameters" on page 4-4
- "Centrifugation" on page 4-10
- "Temperature adaptation during standstill" on page 4-11
- "Short-term centrifugation" on page 4-11
- "Removing the rotor" on page 4-11
- "Turning off the centrifuge" on page 4-12

4

Switch on centrifuge

- 1. Turn on the power switch on the back of the device. The device performs a self-check of its software.
 - a. When the centrifuge lid is closed the following display shows:



The top line shows the speed and the time. Both lines display **0**. The temperature indicator displays the current temperature inside the rotor chamber. The bottom line shows the values for the last run.

b. When the centrifuge lid is open the following display shows:



The top line shows the message lid open. The bottom line shows the values for the last run.

Door opening

c. Press the 🔜 key.



WARNING Do not reach into the crack between the lid and the housing. The lid is drawn shut automatically.

Use the emergency release only for malfunctions and power failures (see "Mechanical emergency lid release" on page 7-2).

Close door

2. Close the door by pressing down on it lightly in the middle or on both sides of it. One lock closes the lid completely.

Note The door should audibly click into place.

Rotor installation

The approved rotors for the Sorvall Legend X1R are listed in section "Rotor selection" on page 1-6. Use only the rotors and accessories from this list in the centrifuge.



CAUTION Unapproved or incorrectly combined accessories can cause serious damage to the centrifuge.

The centrifuge is equipped with a Thermo ScientificTM Auto-LockTM locking system. This system is used to automatically lock the rotor to the centrifuge spindle. The rotor does not have to be bolted on to the centrifuge spindle.

Proceed as follows:

1. Open the lid of the centrifuge and if necessary remove any dust, foreign objects or residue from the chamber.

Auto-Lock and o-ring must be clean and undamaged.

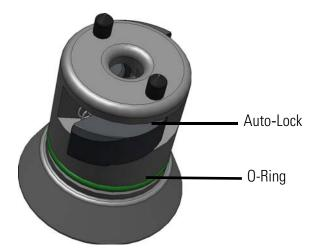


Figure 4-1. Auto-Lock

2. Place the rotor over the centrifuge spindle and let it slide slowly down the centrifuge spindle. The rotor clicks automatically into place.



CAUTION Do not force the rotor onto the centrifuge spindle. If the rotor is very light, then it may be necessary to press it onto the centrifuge spindle with a bit of pressure.

3. Check if the rotor is properly installed by lifting it slightly on the handle. If the rotor can be pulled up, then it must be reclamped to the centrifuge spindle.



WARNING If the rotor cannot be properly locked in place after several attempts, then the Auto-Lock is defective and you are not permitted to operate the rotor. Check for any damage to the rotor: Damaged rotors must not be used. Keep the centrifuge spindle area of the rotor clear of objects.



CAUTION Check that the rotor is properly locked on the centrifuge spindle before each use by pulling it at its handle.

4. If available close the rotor with the rotor lid.



See the information in the rotor instruction manual.

Be sure to check all sealings before starting any aeorsol-tight applications.

5. The display shows:

Entering parameters

Acceleration curve

The Sorvall Legend X1/X1R offers you a total of 9 acceleration and 10 braking curves with which samples and gradients can be centrifuged.

After the centrifuge is turned on, the last running profile selected is shown.

1. Press the <u>selection</u> key once in order to open the acceleration profile selection menu. The display shows the message **Set acceleration**.



- 2. Select the desired profile by entering a digit between 1-9 on the numeric pad. Curve number 1 is the slowest and curve number 9 the fastest.
- 3. Confirm your entry by pressing errer

Braking curve

 Press the kay twice in order to open the braking curve selection menu. The display shows the message Set deceleration:



- 2. Select the desired curve by entering a digit between 1-9 on the numeric pad. Curve number 1 is the slowest and curve number 10 the fastest.
- 3. Confirm your entry by pressing

Preselecting Speed / RCF

1. Press the setto key.

The display shows the **RPM** or the "RCF"-value depending on the display setting. Press the key to toggle between the two modes.



- 2. Enter the desired value using the numeric pad. The digits show in sequential order.
- Confirm your entry by pressing
 Your entry will be automatically confirmed if you do not press any key for 5 seconds.

Note If an extremely low RCF value has been selected, it will be corrected automatically if the resulting speed is less than 300 rpm. This is because 300 rpm is the lowest selectable speed.

Explanation of RCF value

The relative centrifugal force (RCF) is given as a multiple of the force of gravity g. It is a unitless numerical value which is used to compare the separation or sedimentation capacity of various devices, since it is independent of the type of device. Only the centrifuging radius and the speed come into play in it:

$$\mathsf{RCF} = 11, 18 \times \left\langle \frac{\mathsf{n}}{1000} \right\rangle^2 \times \mathsf{r}$$

r = centrifuging radius in cm

n = rotational speed in rpm

The maximum RCF value is related to the maximum radius of the tube opening.

Remember that this value is reduced depending on the tubes and adapters used.

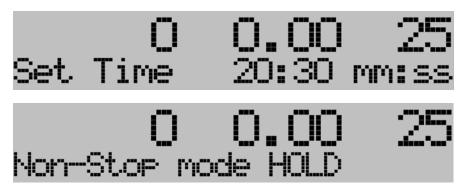
This can be accounted for in the calculation above if required.

Run time preselection

 Press the me key in order to open the runtime selection menu. The display shows the message Set Time. Enter the desired runtime in hh.mm



2. Press the we repeatedly to switch over to mm:ss or Hold mode respectively.



- 3. Enter the desired runtime using the numeric pad. The digits show in sequential order.
- Confirm your entry by pressing
 Your entry will be automatically confirmed if you do not press any key for 5 seconds.

While the centrifuge is in operation, two different values can be displayed. The total run time and the run time after reaching the maximum speed.

Press the 💿 key to toggle between the two display modes.

Continuous operation

During continuous operation, the centrifuge will continue running until you stop it manually with the key.

Run time mode

Use the 💿 key to set the run time counter start.

AT START

When AT START is lit, the run time will be counted right from pressing the key.

AT SPEED

When AT SPEED is lit, the run time will be counted once the preset speed or the preset RCF-value respectively have been reached. In this case, the selected acceleration curve has no effect on the run time.

Preselecting the temperature

You can preselect temperatures between -10 °C and +40 °C.

While the centrifuge is in operation, two different values can be displayed. The sample temperature and the chamber temperature. Press the skey to toggle between the two display modes.

AIR

When AIR is lit, the temperature inside the rotor chamber will be controlled. This setting is recommended when using the centrifuges pretemp feature.

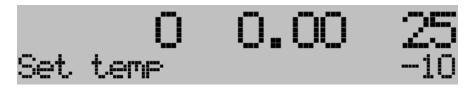
SAMPLE

When SAMPLE is lit, the temperature of the sample will be controlled.

To set the temperature, proceed as follows:

1. Press the key *me* in order to open the temperature selection menu.

The display shows the messages **Temperature**. Depending on the display mode setting, the value shown corresponds either to the sample temperature or to the inside chamber temperature. Press the state to toggle between the two modes.



- 2. Enter the desired temperature using the numeric pad. The digits show in sequential order.
- Confirm your entry by pressing one.
 Your entry will be automatically confirmed if you do not press any key for 5 seconds.

Prewarming or precooling the centrifuge

For setting the pretemp value for the centrifuge proceed as follows:

1. Hold down the wey for a minimum of three seconds to open the temperature selection menu. The display shows the message **Set PreTemp**.



- 2. Enter the desired temperature using the numeric pad. The digits show in sequential order.
- 3. Confirm your entry by pressing area. Your entry will be automatically confirmed if you do not press any key for 5 seconds.



4. The display shows:



5. Press the key.

The rotor chamber is cooled down or heated up to the preset temperature.



6. Press the 🔜 key.

The display shows the current temperature.



Bucket selection

Bucket selection is only possible for swing-out rotors. The bucket code corresponds to the last four digits of the bucket catalog number.

1. Press the 🔳 key.

The display shows the following message:



2. Press the skey repeatedly until the bucket being used is displayed. Your entry will be automatically confirmed if you do not press any key for 5 seconds.

Bucket radius

Having entered the bucket code, you may want to specify the radius.

When you use adapters, you will have to specify the radius. For appropriate values please refer to the instruction manual for the rotor.

1. Press to confirm your selection and open the radius input menu.



The default value is marked accordingly. You can either confirm the default value by pressing or enter a different radius.

2. Enter the radius using the numeric pad. The digits show in sequential order. 3. Press to confirm the new radius

Setting up a program

- 1. Enter the program parameters.
- 2. As described above, please enter braking and acceleration profile, speed/ RCF value, run time, temperatures, bucket code and bucket radius.
- 3. Press any of the program store keys for 4 seconds.

Configuring, storing and protecting a program

Programs can be stored and retrieved later by means of the assigned number.

- 1. Enter the program parameters as described above.
- Press the program memory key to load any of the programs 1-5.
 Press the key in order to load any of the programs 1-99 by means of the numeric pad.
 Enter the program number and confirm with .
- 3. You may want to enter a passcode to protect the data. After four seconds the display shows the message **Program saved** or **Program locked**.



Overwriting a passcode-protected program

If you wish to overwrite a passcode-protected program proceed as follows:

- 1. Select the program you wish to change.
- Press the program memory key to load any of the programs 1-5.
 Press the key in order to load any of the programs 1-99 by means of the numeric pad. Enter the program number and confirm with .



3. Enter the passcode. If you enter the wrong password, the following message appears:



The display restores the current values. Return to step 1.

Centrifugation

Once the rotor has been properly installed, the main switch has been turned on and the lid has been closed, you can start centrifuging.

Starting centrifuge program

Press the key on the control panel. The centrifuge accelerates to the preset speed with the time display active.

If the speed setting is higher than the maximum permissible speed or RCF-value for the particular rotor, then the display will show the message **CHECK SET SPEED**. The centrifuge decelerates. Please enter a new value.

You cannot open the lid as long as the centrifuge is running.

Imbalance indicator

If a load is imbalanced, this will be indicated at speed higher than approx. 300 rpm by the message **Imbalance load**.

The run will terminate.

Check the loading and start the centrifuge once again. See the information on proper loading in the rotor instruction manual. For information on troubleshooting, see section "Troubleshooting by user" on page 7-3.

Stopping the centrifugation program

With preset run time

Usually the run time is preset and you only have to wait until the centrifuge stops automatically when the preset time limit expires.

As soon as the speed drops to zero, the message **End** will appear in the display. By pressing the **a** key, you can open the door and remove the centrifuge material.

You can also stop the centrifuging program manually at any time by pressing the 🔜 key.

Continuous operation

If you selected continuous operation (see "Continuous operation" on page 4-6), you will have to stop the centrifuge manually. Press the set was not been control panel. The centrifuge will be decelerated at the designated rate. The message End will illuminate, and after pressing the set, the lid will open and you can remove the centrifuged material.

Temperature adaptation during standstill

The temperature cannot be adapted until the rotor has been positively identified. For this, the centrifuge must have accelerated to over 500 rpm. The speed display will then show **End**.

When the rotor is not recognized (lid closed and key not yet pressed, speed display **0**), the centrifuge responds by ensuring that the sample cannot freeze regardless of the rotor being used.

Short-term centrifugation

For short-term centrifuging, the Sorvall Legend X1R has a PULSE- function.

By holding down the 🚵 key, spinning will start and continue until the key is let go.

The centrifuge accelerates and brakes at maximum power. Any rpm or RCF entered beforehand is overridden.

Note The centrifuge accelerates to maximum speed, regardless of which rotor was installed.

Check carefully whether you have to maintain a certain speed for your application.

During the acceleration process, time is counted forwards in seconds. The reading stays displayed until the centrifuge lid is opened.

Removing the rotor

To remove the rotor, proceed as follows:

- 1. Open the centrifuge lid.
- 2. Grab the rotor handle with both hands and press against the green Auto-Lock key. At the same time, pull the rotor directly upwards with both hands and remove it from the centrifuge spindle. Make sure not to tilt the rotor while doing this.



Figure 4-2. Press Auto-Lock key

Aerosol-tight rotors

When using an aerosol tight lid the rotor can only be removed with the lid closed. This is to protect you and the samples.



CAUTION Rotors supplied with a lid for aerosol-tight applications come with a mandrel, which belongs to the Auto-Lock. Be sure not to place the lid onto this mandrel to prevent it from being damaged.

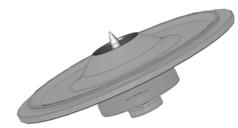


Figure 4-3. Auto-Lock-lid for aerosol-tight rotors

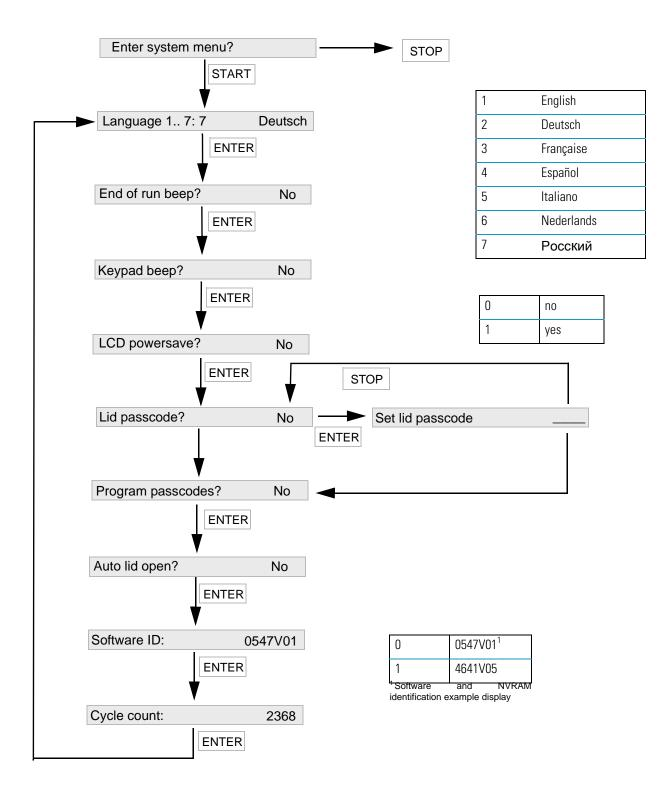


WARNING Mind the Auto-Lock-mandrel inside the lid. Do not touch the mandrel.

Turning off the centrifuge

1. To turn off the centrifuge put the mains switch to "0".

System menu



Description

Enter system menu

To enter the system menu hold down any of the keys when switching the centrifuge on.

Press the me key in order to navigate through the system menu.

Press the 🔜 key to quit the system menu.

Language

In order to change the language in the display, press the number next to the desired language in the diagram on the numeric pad.

Press key to confirm.

End of run beep

Press 1 (yes) on the numeric pad if the centrifuge should make beep after the run. Otherwise press 0 (no).

Press key to confirm.

Keypad beep

Press 1 (yes) on the numeric pad if the centrifuge should make beep when pressing any key. Otherwise press 0 (no).

Press key to confirm.

LCD powersave

Press 1 (yes) on the numeric pad if the centrifuge should enter a powdersave mode after the run. Otherwise press 0 (no).

Press key to confirm.

Lid passcode

Press 1 (yes) on the numeric pad if the centrifuge door should be protected with a passcode. Enter passcode.

If the door should not be protected with a passcode press 2 (no) on the numeric pad.

Press key to confirm.

Program passcodes

Press 1 (yes) on the numeric pad if the centrifuge programs should be protected with a passcode. Otherwise press 0 (no).

Press key to confirm.

Auto lid open

Press 1 (yes) on the numeric pad if the centrifuge should open after the run. Otherwise press 0 (no).

Press key to confirm.

Software ID

Here you fine the current software version.

Press were key to confirm.

Cycle counter

Here you fine the current numbers of cycles.

Press were key to confirm.

Maintenance and care

Contents

- "Cleaning intervals" on page 6-2
- "Cleaning" on page 6-2
- "Disinfection" on page 6-4
- "Decontamination" on page 6-5
- "Autoclaving" on page 6-5
- "Service of Thermo Fisher Scientific" on page 6-6

6



CAUTION During maintenance operations, always disconnect the centrifuge from the power.

Cleaning intervals

For the sake of personal, environmental, and material protection, it is your duty to clean and if necessary disinfect the centrifuge on a regular basis.

Maintenance	Recommended interval
Cabinet / Housing	every month or when polluted
Rotor chamber	daily or when polluted
Rotor	daily or when polluted
Accessories	daily or when polluted
Cabinet	once per month
Ventilation holes	every six months by a service technician



CAUTION Refrain from using any other cleaning or decontamination procedure than those recommended here, if you are not entirely sure that the intended procedure is safe for the equipment.

Use only approved cleansers.

If in doubt, contact Thermo Fisher Scientific.

Cleaning

When cleaning centrifuge mind the following:

- Use warm water with a neutral solvent.
- Never use caustic cleaning agents such as soap suds, phosphoric acid, bleaching solutions or scrubbing powder.
- Rinse the cavities out thoroughly.
- Use a soft brush without metal bristles to remove stubborn residue.
- Afterwards rinse with distilled water.
- Place the rotors on a plastic grate with their cavities pointing down.
- If drying boxes are used, the temperature must never exceed 50 °C, since higher temperatures could damage the material and shorten the lifetime of the parts.
- Use only disinfectants with a pH of 6-8.
- Dry aluminum parts off with a soft cloth.

- After cleaning, treat the entire surface of aluminum parts with corrosion protection oil (7000 9824). Also treat the cavities with oil.
- Store the aluminum parts at room temperature or in a cold-storage room with the cavities pointing down.



CAUTION Before using any cleaning or decontamination methods except those recommended by the manufacturer, users should check with the manufacturer that the proposed method will not damage the equipment.

Clean centrifuge and accessories as follows:

- 1. Open the centrifuge.
- 2. Turn off the centrifuge.
- 3. Pull out the power supply plug.
- 4. Grasp the rotor with both hands and lift it vertically off the centrifuge spindle.
- 5. Remove the centrifuge tubes and adapters.
- 6. Use a neutral cleaning agent with a pH value between 6 and 8 for cleaning.
- 7. Use a small nylon brush to clean the cavities of fixed angle rotors and buckets.



CAUTION Do not use metal wire brushes or other tools that may scratch the anodized surface of the buckets or rotors.

- 8. Dry all of the rotors and accessories after cleaning with a cloth or in a warm air cabinet at a maximum temperature of 50 °C.
- After cleaning, treat the entire surface of aluminum parts with corrosion protection oil (7000 9824). Also treat the cavities with oil.
- Tread the bold of the swing out rotor with bold grease (75003786).



CAUTION When cleaning, do not allow liquids, especially organic solvents, to get on the drive shaft or the bearings of the centrifuge.

Organic solvents break down the grease in the motor bearing. The drive shaft could freeze up.

After some applications their might be ice in the rotor chamber. Let the ice melt and drain it off. Clean the rotor chamber as described above.

Disinfection

Disinfect the centrifuge immediately whenever infectious material has spilled during centrifugation.



WARNING Infectious material can get into the centrifuge when a tube breaks or as a result of spills. Keep in mind the risk of infection when touching the rotor and take all necessary precautions.

In case of contamination, make sure that others are not put at risk. Decontaminate the affected parts immediately.

Take other precautions if need be.

The rotor chamber and the rotor should be treated preferably with a neutral disinfectant.



CAUTION Before using any cleaning or decontamination methods except those recommended by the manufacturer, users should check with the manufacturer that the proposed method will not damage the equipment.

Observe the safety precautions and handling instructions for the cleaning agents used.

Contact the Service Department of Thermo Fisher Scientific for questions regarding the use of other disinfectants.

Disinfect the rotor and accessories as follows:

- 1. Open the centrifuge.
- 2. Turn off the centrifuge.
- 3. Pull out the power supply plug.
- 4. Grasp the rotor with both hands and lift it vertically off the centrifuge spindle.
- 5. Remove the centrifuge tubes and adapters and dispose of them or disinfect them.
- 6. Treat the rotor and accessories according to the instructions for the disinfectant. Adhere strictly to the given application times.
- 7. Be sure the disinfectant can drain off the rotor.
- 8. Rinse the rotor and accessories thoroughly with water.
- 9. Dispose of the disinfectant according to the applicable guidelines.
- 10. Dry all of the rotors and accessories after cleaning with a cloth or in a warm air cabinet at a maximum temperature of 50 $^{\circ}$ C.
 - After cleaning, treat the entire surface of aluminum parts with corrosion protection oil (7000 9824). Also treat the cavities with oil.
 - Tread the bold of the swing out rotor with bold grease (75003786).

Decontamination

Decantaminate the centrifuge immediately whenever radioactive material has spilled during centrifugation.



WARNING Radioactive material can get into the centrifuge when a tube breaks or as a result of spills. Keep in mind the risk of infection when touching the rotor and take all necessary precautions.

In case of contamination, make sure that others are not put at risk. Decontaminate the affected parts immediately. Take other precautions if need be.



CAUTION Before using any cleaning or decontamination methods except those recommended by the manufacturer, users should check with the manufacturer that the proposed method will not damage the equipment.

For general radioactive decontamination use a solution of equal parts of 70% ethanol, 10% SDS and water.

- 1. Open the centrifuge.
- 2. Turn off the centrifuge.
- 3. Pull out the power supply plug.
- 4. Grasp the rotor with both hands and lift it vertically off the centrifuge spindle.
- 5. Remove the centrifuge tubes and adapters and dispose of them or disinfect them.
- 6. Rinse the rotor first with ethanol and then with de-ionized water.
- Adhere strictly to the given application times.
- 7. Be sure the decontamination solution can drain off the rotor.
- 8. Rinse the rotor and accessories thoroughly with water.
- 9. Dispose of the decontamination solution according to the applicable guidelines.
- 10. Dry all of the rotors and accessories after cleaning with a cloth or in a warm air cabinet at a maximum temperature of 50 $^{\circ}$ C.
 - After cleaning, treat the entire surface of aluminum parts with corrosion protection oil (7000 9824). Also treat the cavities with oil.
 - Tread the bold of the swing out rotor with bold grease (75003786).

Autoclaving

- 1. Before autoclaving clean rotor and accessories and described above.
- 2. Place the rotor on a flat surface.
- Rotors and adapter can be autoclaved at 121 °C.

• The maximum permissible autoclave cycle is 20 minutes at 121 °C.

Note No chemical additives are permitted in the steam.



CAUTION Never exceed the permitted temperature and duration when autoclaving. If the rotor shows signs of corrosion or wear, it must be replaced.

Service of Thermo Fisher Scientific

Thermo Fisher Scientific recommends having the centrifuge and accessories serviced once a year by an authorized service technician. The service technician check the following:

- the electrical equipment
- the suitability of the set-up site
- the lid lock and the safety system
- the rotor

technician.

• the fixation of the rotor and the drive shaft

Thermo Fisher Scientific offers inspection and service contracts for this work. Any necessary repairs are performed for free during the warranty period and afterwards for a charge. This is only valid if the centrifuge has only been maintained by a Thermo Fisher Scientific service

Sorvall Legend X1R

6-6

Troubleshooting

Contents

- "Mechanical emergency lid release" on page 7-2
- "Troubleshooting by user" on page 7-3
- "When to contact customer service" on page 7-6

Mechanical emergency lid release

During a power failure, you will not be able to open the centrifuge lid with the regular electric lid release. A mechanical override is provided to allow sample recovery in the case of an emergency. However, this should be used only in emergencies and after the rotor has come to a complete stop.



WARNING The rotor can still be spinning at high speed. If touched, it can cause serious injuries.

Always wait a few minutes until the rotor has come to a stop without braking. The brake does not work when there is no current. The braking process lasts much longer than usual.

Proceed as follows:

1. Make sure the rotor has stopped (view port in the lid).



WARNING Never use your hand or other tools to brake the rotor.

- 2. Pull out the power supply plug.
- 3. On the left side of the housing is one white plastic plugs which you can pry out of the side plate with a screwdriver or a knife.
 Pull the release cord attached to it to trigger the mechanical lid release. The lid will open and the

Pull the release cord attached to it to trigger the mechanical lid release. The lid will open and the samples can be removed.

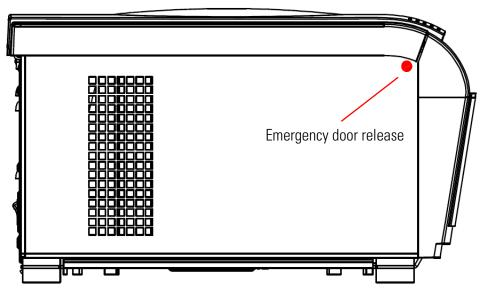


Figure 7-1. Emergency door release

4. Push the cord back into the centrifuge and mount the plug.

Reconnect the centrifuge once the power has been restored. Switch on the centrifuge. Press the 🔜 key to have the door locks operative again.

Troubleshooting by user



If problems occur other than those listed in this table, the authorized customer service representative must be contacted.

Failure message	Problem with centrifuge	Possible causes and cures
Chamber overtemp	The centrifuge cannot be operated. The run does not start or the centrifuge runs down without being braked.	Overheating in chamber. Check the function of the refrigeration unit. Clean the air inlet for the condenser.
	Ū	Restart the centrifuge. If an error message appears again, inform your service technician.
Incorrect bucket ID	The centrifuge cannot be operated. The run does not start or the centrifuge runs down without being braked.	Bucket code is undefined for the rotor detected, check the set points for the given bucket code. Is it permitted to use the current bucket in the rotor currently mounted?
		Restart the centrifuge. If an error message appears again, inform your service technician.
Unapproved rotor	The centrifuge cannot be operated.	Rotor code is not in the rotor table.
	The run does not start or the centrifuge runs	Is it permitted to use the rotor currently mounted in this device?
	down without being braked.	Restart the centrifuge. If an error message appears again, inform your service technician.
Rotor ID failure	The centrifuge cannot be operated.	The rotor could not be identified.
	The run does not start or the centrifuge runs down without being braked.	Check to see if the rotor is properly installed. Restart the centrifuge. If an error message appears again, inform your service technician.
Lid blocked	Centrifuge does not open	Restart the centrifuge.
		The emergency lid release enables you to retrieve your samples
		If an error message appears again, inform your service technician.
Motor overtemp	The centrifuge cannot be operated.	Restart the centrifuge.
	The run does not start or the centrifuge runs down without being braked.	If an error message appears again, inform your service technician.
PCB overtemp	The centrifuge cannot be operated.	Restart the centrifuge.
	The run does not start or the centrifuge runs down without being braked.	If an error message appears again, inform your service technician.

Failure message	Problem with centrifuge	Possible causes and cures
Error lid lock	The centrifuge cannot be operated.	The lid opens while the device is running.
	The run does not start or the centrifuge runs	Close the lid and restart the centrifuge.
	down without being braked.	If an error message appears again, inform your service technician.
Imbalanced load	The centrifuge cannot be operated.	Imbalance detected.
	The run does not start or the centrifuge runs	Check the load placed in the rotor.
	down without being braked.	Check that the rotor cross bolts are well greased.
		Restart the centrifuge.
		If an error message appears again, inform your service technician.
Check Set Speed	The centrifuge cannot be operated.	The set point speed is higher than the maximum rotor speed.
	The run does not start or the centrifuge runs	Correct the value.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-01 - E-12	The centrifuge cannot be operated.	Error during the self-test of the centrifuge program and the
	The run does not start or the centrifuge runs	electronics.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-13	The centrifuge cannot be operated.	The check sum in the data memory is incorrect.
	The run does not start or the centrifuge runs	The software corrects errors automatically.
	down without being braked.	Check the values of the set point settings, etc.
		Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-15-E-16	Temperature sensor broken / controller	Malfunction in the temperature detection.
	defective	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-17	Speed for rotor detection exceeded	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-21 - E-22	The centrifuge cannot be operated.	The rotor could not be identified.
	The run does not start or the centrifuge runs	Check to see if the rotor is properly installed.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-23	The centrifuge cannot be operated.	The speed control measurement returned a different result.
	The run does not start or the centrifuge runs	Restart the centrifuge.
	down without being braked.	If an error message appears again, inform your service technician.

Failure message	Problem with centrifuge	Possible causes and cures
E-25-E-27	The centrifuge cannot be operated.	Time has expired for the lid lock drive while opening the lid.
	The run does not start or the centrifuge runs	Restart the centrifuge.
	down without being braked.	If an error message appears again, inform your service technician.
E-28	The centrifuge cannot be operated.	Error during the self-test of the centrifuge program and the
	The run does not start or the centrifuge runs	electronics.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-29	The centrifuge cannot be operated.	Check wether you selected the right bucket.
	The run does not start.	Is it easy to turn the rotor when the lid is open? Does the rotor rub against the device?
		Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-30	The centrifuge cannot be operated.	Restart the centrifuge.
	The run does not start or the centrifuge runs down without being braked.	If an error message appears again, inform your service technician.
E-33	The centrifuge cannot be operated.	Overpressure in the refrigeration unit.
	The run does not start or the centrifuge runs	Clean the air inlet for the condenser.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-34-E-36	The centrifuge cannot be operated.	Error during the self-test of the centrifuge program and the
	The run does not start or the centrifuge runs	electronics.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-40	The centrifuge cannot be operated.	The centrifuge accelerates too slowly.
	The run does not start or the centrifuge runs down without being braked.	Is the rotor properly installed? Check wether you selected the right bucket. Is it easy to turn the rotor when the lid is open? Does the rotor rub against the device?
		Restart the centrifuge.
		If an error message appears again, inform your service technician.
E-41-E-74	The centrifuge cannot be operated. The run does not start or the centrifuge runs	Error during the self-test of the centrifuge program and the electronics.
	down without being braked.	Restart the centrifuge.
		If an error message appears again, inform your service technician.

When to contact customer service

If you need to contact customer service, please provide the order no. and the serial no. of your device. This information can be found on the back near the inlet for the power supply cable.

To identify the software version, proceed as follows:

- 1. Hold down any of the keys and then switch on the centrifuge. You enter the system menu.
- 2. Press the 📩 key.
- 3. Press and hold the event key, until the following message is displayed:



4. Communicate the software version to the service technician.

Chemical compatibility chart

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING For ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL	NATON	PET*, POLYCLEAR, CLEARCRIMP	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYTHERMIDE	POLYRTHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A, TEFLON	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON	VITON
2-mercaptoethanol		S	S	U	-	S	Μ	S	-	S	U	S	S	U	S	S	-	S	S	S	S	U	S	S	S	S	S	S
Acetaldehyde		S	-	U	U	-	-	-	Μ	-	U	-	-	-	Μ	U	U	U	Μ	Μ	-	Μ	S	U	-	S	-	U
Acetone		Μ	S	U	U	S	U	Μ	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	Μ	Μ	S	U	U
Acetonitrile		S	S	U	-	S	Μ	S	-	S	S	U	S	U	Μ	U	U	-	S	Μ	U	U	S	S	S	S	U	U
Alconox		U	U	S	-	S	S	S	-	S	S	S	S	S	S	Μ	S	S	S	S	S	S	S	S	S	S	S	U
Allyl Alcohol		-	-	-	U	-	-	S	-	-	-	-	S	-	S	S	Μ	S	S	S	-	Μ	S	-	-	S	-	-
Aluminum Chloride		U	U	S	S	S	S	U	S	S	S	S	Μ	S	S	S	S	-	S	S	S	S	S	Μ	U	U	S	S
Formic Acid (100%)		-	S	Μ	U	-	-	U	-	-	-	-	U	-	S	Μ	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		Μ	S	U	S	S	S	S	S	S	S	S	S	S	S	U	U	-	S	S	S	S	S	S	Μ	S	S	S
Ammonium Hydroxide (10%)		U	U	S	U	S	S	Μ	S	S	S	S	S	-	S	U	Μ	S	S	S	S	S	S	S	S	S	Μ	S
Ammonium Hydroxide (28%)		U	U	S	U	S	U	Μ	S	S	S	S	S	U	S	U	Μ	S	S	S	S	S	S	S	S	S	Μ	S
Ammonium Hydroxide (conc.)		U	U	U	U	S	U	Μ	S	-	S	-	S	U	S	U	U	S	S	S	-	Μ	S	S	S	S	-	U
Ammonium Phosphate		U	-	S	-	S	S	S	S	S	S	S	S	-	S	S	Μ	-	S	S	S	S	S	S	Μ	S	S	S
Ammonium Sulfate		U	Μ	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	-	Μ	U	-	-	S	S	-	Μ	-	S	-	Μ	S	S	S	S	М	-	-	-	U	-	S	-	Μ
Aniline		S	S	U	U	S	U	S	Μ	S	U	U	U	U	U	U	U	-	S	М	U	U	S	S	S	S	U	S
Sodium Hydroxide (<1%)		U	-	Μ	S	S	S	-	-	S	Μ	S	S	-	S	Μ	Μ	S	S	S	S	S	S	Μ	S	S	-	U
Sodium Hydroxide (10%)		U	-	Μ	U	-	-	U	-	Μ	Μ	S	S	U	S	U	U	S	S	S	S	S	S	Μ	S	S	-	U
Barium Salts		Μ	U	S	-	S	S	S	S	S	S	S	S	S	S	S	Μ	-	S	S	S	S	S	S	Μ	S	S	S
Benzene		S	S	U	U	S	U	Μ	U	S	U	U	S	U	U	U	Μ	U	Μ	U	U	U	S	U	U	S	U	S
Benzyl Alcohol		S	-	U	U	-	-	Μ	Μ	-	Μ	-	S	U	U	U	U	U	U	U	-	Μ	S	Μ	-	S	-	S
Boric Acid		U	S	S	Μ	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		М	-	S	-	S	S	S	-	S	S	S	S	-	S	S	-	-	S	S	S	S	S	S	М	S	S	S

A

CHEMICAL	MATERIAL	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL	NYLON	PET*, POLYCLEAR, CLEARCRIMP	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYTHERMIDE	POLYRTHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	rulon a, teflon	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON	VITON
Cesium Bromide	Μ	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Cesium Chloride	М	S	S	U	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Cesium Formate	М	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Cesium lodide	М	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Cesium Sulfate	М	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Chloroform	U	U	U	U	S	S	Μ	U	S	U	U	Μ	U	Μ	U	U	U	Μ	Μ	U	U	S	U	U	U	Μ	S
Chromic Acid (10%)	U	-	U	U	S	U	U	-	S	S	S	U	S	S	Μ	U	Μ	S	S	U	Μ	S	Μ	U	S	S	S
Chromic Acid (50%)	U	-	U	U	-	U	U	-	-	-	S	U	U	S	Μ	U	Μ	S	S	U	Μ	S	-	U	Μ	-	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	Μ	S	Μ	U	Μ	Μ	S	U	Μ	Μ	U	S
Deoxycholate	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	М	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	Μ	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	Μ	U
Diethyl Ketone	S	-	U	U	-	-	Μ	-	S	U	-	S	-	Μ	U	U	U	Μ	Μ	-	U	S	-	-	S	U	U
Diethylpyrocarbonate	S	S	U	-	S	S	S	-	S	S	U	S	U	S	U	-	-	S	S	S	Μ	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	Μ	S	U	U	S	S	Μ	Μ	S	U	U	S	U	Μ	U	U	-	Μ	Μ	Μ	U	S	S	S	S	U	U
Ferric Chloride	U	U	S	-	-	-	Μ	S	-	Μ	-	S	-	S	-	-	-	S	S	-	-	-	Μ	U	S	-	S
Acetic Acid (Glacial)	S	S	U	U	S	S	U	Μ	S	U	S	U	U	U	U	U	Μ	S	U	Μ	U	S	U	U	S	-	U
Acetic Acid (5%)	S	S	Μ	S	S	S	Μ	S	S	S	S	S	Μ	S	S	S	S	S	S	S	Μ	S	S	Μ	S	S	Μ
Acetic Acid (60%)	S	S	U	U	S	S	U	-	S	Μ	S	U	U	Μ	U	S	Μ	S	Μ	S	Μ	S	Μ	U	S	Μ	U
Ethyl Acetate	Μ	Μ	U	U	S	S	Μ	Μ	S	S	U	S	U	Μ	U	U	-	S	S	U	U	S	Μ	Μ	S	U	U
Ethyl Alcohol (50%)	S	S	S	S	S	S	Μ	S	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	Μ	S	Μ	U
Ethyl Alcohol (95%)	S	S	S	U	S	S	Μ	S	S	S	S	S	U	S	U	-	S	S	S	Μ	S	S	S	U	S	Μ	U
Ethylene Dichloride	S	-	U	U	-	-	S	Μ	-	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	Μ	S	Μ	S
Ethylene Oxide Vapor	S	-	U	-	-	U	-	-	S	U	-	S	-	S	Μ	-	-	S	S	S	U	S	U	S	S	S	U
Ficoll-Hypaque	М	S	S	-	S	S	S	-	S	S	S	S	-	S	S	-	S	S	S	S	S	S	S	Μ	S	S	S
Hydrofluoric Acid (10%)	U	U	U	Μ	-	-	U	-	-	U	U	S	-	S	Μ	U	S	S	S	S	Μ	S	U	U	U	-	-
Hydrofluoric Acid (50%)	U	U	U	U	-	-	U	-	-	U	U	U	U	S	U	U	U	S	S	Μ	Μ	S	U	U	U	-	Μ
Hydrochloric Acid (conc.)	U	U	U	U	-	U	U	Μ	-	U	Μ	U	U	Μ	U	U	U	-	S	-	U	S	U	U	U	-	-

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL	Nylon	PET*, POLYCLEAR, CLEARCRIMP	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYTHERMIDE	POLYRTHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A, TEFLON	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON	VITON
Formaldehyde (40%)		Μ	Μ	Μ	S	S	S	S	Μ	S	S	S	S	Μ	S	S	S	U	S	S	Μ	S	S	S	Μ	S	Μ	U
Glutaraldehyde		S	S	S	S	-	-	S	-	S	S	S	S	S	S	S	-	-	S	S	S	-	-	S	S	S	-	-
Glycerol		Μ	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S
Guanidine 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	Μ	U	S	S	U	S	S	Μ	S	U	S	S	U	S
Isobutyl Alcohol		-	-	Μ	U	-	-	S	S	-	U	-	S	U	S	S	Μ	S	S	S	-	S	S	S	-	S	-	S
Isopropyl Alcohol		Μ	Μ	М	U	S	S	S	S	S	U	S	S	U	S	U	Μ	S	S	S	S	S	S	S	Μ	Μ	Μ	S
Iodoacetic Acid		S	S	М	-	S	S	S	-	S	Μ	S	S	Μ	S	S	-	Μ	S	S	S	S	S	Μ	S	S	Μ	М
Potassium Bromide		U	S	S	-	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	Μ	S	S	S
Potassium Carbonate		М	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	Μ	-	S	S	S	S	-	S	U	S	S	S	S	S	S	S	М	U	Μ	S	U
Potassium Hydroxide (conc.)		U	U	М	U	-	-	Μ	-	Μ	S	S	-	U	Μ	U	U	U	S	Μ	-	Μ	U	-	U	U	-	U
Potassium Permanganate		S	S	S	-	S	S	S	-	S	S	S	U	S	S	S	Μ	-	S	Μ	S	U	S	S	Μ	S	U	S
Calcium Chloride		Μ	U	S	S	S	S	S	S	S	S	S	S	S	S	Μ	S	-	S	S	S	S	S	S	Μ	S	S	S
Calcium Hypochlorite		Μ	-	U	-	S	Μ	Μ	S	-	М	-	S	-	S	Μ	S	-	S	S	S	Μ	S	Μ	U	S	-	S
Kerosene		S	S	S	-	S	S	S	U	S	М	U	S	U	Μ	Μ	S	-	Μ	Μ	Μ	S	S	U	S	S	U	S
Sodium Chloride (10%)		S	-	S	S	S	S	S	S	-	-	-	S	S	S	S	S	-	S	S	S	S	-	S	S	Μ	-	S
Sodium Chloride (sat'd)		U	-	S	U	S	S	S	-	-	-	-	S	S	S	S	S	-	S	S	-	S	-	S	S	Μ	-	S
Carbon Tetrachloride		U	U	Μ	S	S	U	Μ	U	S	U	U	S	U	Μ	U	S	S	Μ	Μ	S	Μ	Μ	Μ	Μ	U	S	S
Aqua Regia		U	-	U	U	-	-	U	-	-	-	-	-	U	U	U	U	U	U	U	-	-	-	-	-	S	-	Μ
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		Μ	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	Μ	S	S	S
Mercaptoacetic Acid		U	S	U	-	S	Μ	S	-	S	Μ	S	U	U	U	U	-	S	U	U	S	Μ	S	U	S	S	S	S
Methyl Alcohol		S	S	S	U	S	S	Μ	S	S	S	S	S	U	S	U	Μ	S	S	S	S	S	S	S	Μ	S	Μ	U
Methylene Chloride		U	U	U	U	Μ	S	S	U	S	U	U	S	U	U	U	U	U	Μ	U	U	U	S	S	Μ	U	S	U
Methyl Ethyl Ketone		S	S	U	U	S	S	Μ	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	S	S	S	U	U
Metrizamide		Μ	S	S	-	S	S	S	-	S	S	S	S	-	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Lactic Acid (100%)		-	-	S	-	-	-	-	-	-	М	S	U	-	S	S	S	Μ	S	S	-	Μ	S	Μ	S	S	-	S
Lactic Acid (20%)		-	-	S	S	-	-	-	-	-	М	S	Μ	-	S	S	S	S	S	S	S	Μ	S	Μ	S	S	-	S
N-Butyl Alcohol		S	-	S	U	-	-	S	-	-	S	М	-	U	S	Μ	S	S	S	S	М	М	S	М	-	S	-	S

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING For ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL	NOTAN	PET*, POLYCLEAR, CLEARCRIMP	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYTHERMIDE	POLYRTHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A, TEFLON	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON	VITON
N-Butyl Phthalate		S	S	U	-	S	S	S	-	S	U	U	S	U	U	U	Μ	-	U	U	S	U	S	Μ	Μ	S	U	S
N, N-Dimethylformamide		S	S	S	U	S	Μ	S	-	S	S	U	S	U	S	U	U	-	S	S	U	U	S	Μ	S	S	S	U
Sodium Borate		Μ	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	-	S	S	S	S	S	S	Μ	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	Μ	S	S	S
Sodium Carbonate (2%)		Μ	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	Μ	S	S	Μ	U	S	S	Μ	S	S	S	Μ	S	S	S	S	Μ	S	S	S	Μ	U	S	М	S
Sodium lodide		Μ	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	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	S	Μ	S	S	S
Sodium Sulfide		S	-	S	S	-	-	-	S	-	-	-	S	S	S	U	U	-	-	S	-	-	-	S	S	Μ	-	S
Sodium Sulfite		S	S	S	-	S	S	S	S	Μ	S	S	S	S	S	S	Μ	-	S	S	S	S	S	S	S	S	S	S
Nickel Salts		U	S	S	S	S	S	-	S	S	S	-	-	S	S	S	S	-	S	S	S	S	S	S	Μ	S	S	S
Oils (Petroleum)		S	S	S	-	-	-	S	U	S	S	S	S	U	U	Μ	S	Μ	U	U	S	S	S	U	S	S	S	S
Oils (Other)		S	-	S	-	-	-	S	Μ	S	S	S	S	U	S	S	S	S	U	S	S	S	S	-	S	S	Μ	S
Oleic Acid		S	-	U	S	S	S	U	U	S	U	S	S	Μ	S	S	S	S	S	S	S	S	S	Μ	U	S	Μ	Μ
Oxalic Acid		U	U	Μ	S	S	S	U	S	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	U	Μ	S	S
Perchloric Acid (10%)		U	-	U	-	S	U	U	-	S	Μ	Μ	-	-	Μ	U	Μ	S	Μ	Μ	-	Μ	S	U	-	S	-	S
Perchloric Acid (70%)		U	U	U	-	-	U	U	-	S	U	Μ	U	U	Μ	U	U	U	Μ	Μ	U	Μ	S	U	U	S	U	S
Phenol (5%)		U	S	U	-	S	Μ	Μ	-	S	U	Μ	U	U	S	U	Μ	S	Μ	S	U	U	S	U	Μ	Μ	Μ	S
Phenol (50%)		U	S	U	-	S	U	Μ	-	S	U	Μ	U	U	U	U	U	S	U	Μ	U	U	S	U	U	U	Μ	S
Phosphoric Acid (10%)		U	U	Μ	S	S	S	U	S	S	S	S	U	-	S	S	S	S	S	S	S	S	S	U	Μ	U	S	S
Phosphoric Acid (conc.)		U	U	Μ	Μ	-	-	U	S	-	Μ	S	U	U	Μ	Μ	S	S	S	Μ	S	Μ	S	U	Μ	U	-	S
Physiologic Media (Serum, Urine)		Μ	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	Μ	S	S	S	Μ	S	U	S	S	S	U	S	S	S	S	U	S	U	Μ	S	Μ	S
Pyridine (50%)		U	S	U	U	S	U	U	-	U	S	S	U	U	Μ	U	U	-	U	S	Μ	U	S	S	U	U	U	U
Rubidium Bromide		Μ	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Rubidium Chloride		Μ	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	Μ	S	S	S
Sucrose		Μ	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		Μ	S	S	-	S	S	S	-	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	Μ	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	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	Μ	S	S	S	S

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING For ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLY URETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL	NATON	PET*, POLYCLEAR, CLEARCRIMP	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYTHERMIDE	POLYRTHYLENE	POLYPROPYLENE	POLY SULFONE	POLYVINYL CHLORIDE	RULON A, TEFLON	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON	VITON
Nitric Acid (50%)		U	S	U	Μ	S	U	U	-	S	U	S	U	U	Μ	Μ	U	Μ	Μ	Μ	S	S	S	U	S	S	Μ	S
Nitric Acid (95%)		U	-	U	U	-	U	U	-	-	U	U	U	U	Μ	U	U	U	U	Μ	U	U	S	U	S	S	-	S
Hydrochloric Acid (10%)		U	U	Μ	S	S	S	U	-	S	S	S	U	U	S	U	S	S	S	S	S	S	S	S	U	Μ	S	S
Hydrochloric Acid (50%)		U	U	U	U	S	U	U	-	S	Μ	S	U	U	Μ	U	U	S	S	S	S	Μ	S	Μ	U	U	Μ	Μ
Sulfuric Acid (10%)		Μ	U	U	S	S	U	U	-	S	S	Μ	U	S	S	S	S	S	S	S	S	S	S	U	U	U	S	S
Sulfuric Acid (50%)		Μ	U	U	U	S	U	U	-	S	S	Μ	U	U	S	U	U	Μ	S	S	S	S	S	U	U	U	Μ	S
Sulfuric Acid (conc.)		Μ	U	U	U	-	U	U	Μ	-	-	Μ	U	U	S	U	U	U	Μ	S	U	Μ	S	U	U	U	-	S
Stearic Acid		S	-	S	-	-	-	S	Μ	S	S	S	S	-	S	S	S	S	S	S	S	S	S	Μ	Μ	S	S	S
Tetrahydrofuran		S	S	U	U	S	U	U	Μ	S	U	U	S	U	U	U	-	М	U	U	U	U	S	U	S	S	U	U
Toluene		S	S	U	U	S	S	Μ	U	S	U	U	S	U	U	U	S	U	Μ	U	U	U	S	U	S	U	U	Μ
Trichloroacetic Acid		U	U	U	-	S	S	U	Μ	S	U	S	U	U	S	Μ	-	Μ	S	S	U	U	S	U	U	U	Μ	U
Trichloroethane		S	-	U	-	-	-	Μ	U	-	U	-	S	U	U	U	U	U	U	U	U	U	S	U	-	S	-	S
Trichloroethylene		-	-	U	U	-	-	-	U	-	U	-	S	U	U	U	U	U	U	U	U	U	S	U	-	U	-	S
Trisodium Phosphate		-	-	-	S	-	-	Μ	-	-	-	-	-	-	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
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
Urea		S	-	U	S	S	S	S	-	-	-	-	S	S	S	Μ	S	S	S	S	-	S	S	S	Μ	S	-	S
Hydrogen Peroxide (10%)		U	U	Μ	S	S	U	U	-	S	S	S	U	S	S	S	Μ	U	S	S	S	S	S	S	Μ	S	U	S
Hydrogen Peroxide (3%)		S	Μ	S	S	S	-	S	-	S	S	S	S	S	S	S	S	Μ	S	S	S	S	S	S	S	S	S	S
Xylene		S	S	U	S	S	S	Μ	U	S	U	U	U	U	U	U	Μ	U	Μ	U	U	U	S	U	Μ	S	U	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	U	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
Citric Acid (10%)		Μ	S	S	М	S	S	Μ	S	S	S	S	S	S	S	S	S	Μ	S	S	S	S	S	S	S	S	S	S

*Polyethyleneterephthalate

Key

- S Satisfacto y
- 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.

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

Index

A

Acceleration curve	
Accident prevention	
Aerosol-tight rotors	
Aligning the centrifuge	
Autoclaving	

В

Braking curve	4-4
Bucket radius	
Bucket selection	4-8

C

Care	6-1
Centrifugation	
Cleaning	
Cleaning intervals	
Close door	
Contact Information	B-1
Continuous operation	4-6, 4-11
Control panel	
Customer service	

D

Decontamination	6-5
Disinfection	6-4
Door opening	4-2

Ε

Emergency door release	7-2
Entering parameters	
Explanation of RCF value	

F

Failure message7-3
Features1-5

mbalance indicator4-10	

Ι

K

Keys		3-2
------	--	-----

L

```
Location ......2-2
```

Μ

Mains Connection	2-3
Mains supply	1-5, 2-4
Maintenance	6-1

0

Operation	4-1
Overwriting a passcode-protected program	4-9
Overwriting a program	4-9

Ρ

Passcode	
Precautions	v
Preface	iii
Preselecting RCF	4-4
Preselecting speed	4-4
Preselecting the temperature	4-6
Preset run time	4-10
Prewarming or precooling	4-7
Prewarming or precooling the centrifuge .	

R

1
2
6
5
•

S

Safety features	1-2
Safety zone	
Select program	
Short-term centrifugation	
Starting centrifuge program	
Stopping the centrifugation program	
Storage	2-4
Switch on centrifuge	

Т

Technical Data	1-3
Temperature adaptation	
Transporting the centrifuge	
Troubleshooting	

thermo scientific



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

thermofisher.com/centrifuge

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries unless otherwise indicated.

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 Sigma-Aldrich Co. LLC. Valox is a registered trademark of General Electric Co.

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. Pictures of the user interface within the manual are showing the English version as example.

Australia +61 39757 4300 Austria +43 1 801 40 0 Belgium +32 9 272 54 82 China +800 810 5118, +400 650 5118 France +33 2 2803 2180 Germany national toll free 0800 1 536 376 Germany international +49 6184 90 6000 India toll free +1800 22 8374 India +91 22 6716 2200 Italy +39 02 95059 552 Japan +81 3 5826 1616 Korea +82 2 2023 0600 Netherlands +31 76 579 55 55 New Zealand +64 9 980 6700 Nordic/Baltic/CIS countries +358 10 329 2200 Russia +7 812 703 42 15, +7 495 739 76 41 Singapore +82 2 3420 8700 Spain/Portugal +34 93 223 09 18 Switzerland +41 44 454 12 12 UK/Ireland +44 870 609 9203 USA/Canada +1 866 984 3766

Other Asian Countries +852 3107 7600 **Countries not listed** +49 6184 90 6000





