



sartorius

Operating Manual

Centrisart[®] A-14

Laboratory Centrifuge from serial no. 157613



© Copyright by
Sartorius Lab Instruments GmbH & Co. KG
Weender Landstraße 94-108
37075 Göttingen
Germany

Phone: +49.551.308.0
Fax: +49.551.308.3289
www.sartorius.com

1	General information	9
1.1	Importance of the operating manual	9
1.2	Intended use.....	9
1.3	Warranty and liability	9
1.4	Copyright	10
1.5	Standards and regulations	10
1.6	Scope of supply	10
2	Layout and mode of operation.....	11
2.1	Layout of the centrifuge	11
2.1.1	Functional and operating elements	11
2.1.2	Name plate	12
2.2	Mode of operation	13
2.2.1	Centrifugation principle	13
2.2.2	Area of application.....	13
2.2.2.1	Speed, radius, and relative centrifugal force	14
2.2.2.2	Density.....	14
3	Safety	15
3.1	Marking of the unit	15
3.2	Explanation of the symbols and notes	16
3.3	Responsibility of the operator	17
3.4	Operating personnel	17
3.5	Informal safety instructions	17
3.6	Safety instructions	18
3.6.1	Electrical safety	18
3.6.2	Mechanical safety.....	18
3.6.3	Fire prevention	19
3.6.4	Chemical and biological safety.....	19
3.6.5	Safety instructions for centrifugation	20
3.6.6	Resistance of plastics.....	20
3.6.7	Service life of rotors and accessories.....	21
3.7	Safety devices.....	21
3.7.1	Lid lock device	21
3.7.2	System check.....	21
3.8	Measures in the event of hazards and accidents	21
3.9	Remaining hazards.....	22
4	Storage and transport	23
4.1	Dimensions and weight.....	23
4.2	Storage conditions	23
4.3	Notes on transport	23
4.4	Packaging.....	24
4.5	Transport safety device	24

5	Set-up and connection	25
5.1	Installation site.....	25
5.2	Power supply.....	25
5.2.1	Type of connection.....	25
5.2.2	Customer-provided fuses	25
6	Using the centrifuge.....	26
6.1	Initial start-up	26
6.2	Switching the centrifuge on	26
6.2.1	Opening and closing the lid.....	26
6.2.2	Installation of rotors and accessories	27
6.2.2.1	Installation of the rotor	27
6.2.2.2	Installation of accessories.....	28
6.2.2.3	Tubes	28
6.3	Control System "Spincontrol Basic"	29
6.3.1	User interface	29
6.3.2	Display.....	29
6.3.3	Starting a centrifugation run	30
6.3.4	Interrupting a centrifugation run.....	30
6.3.5	Interrupting a deceleration process.....	30
6.3.6	Speed / Relative centrifugal force (RCF).....	30
6.3.6.1	Changing the speed/RCF value during centrifugation.....	30
6.3.7	Runtime	31
6.3.7.1	Changing the runtime during centrifugation.....	31
6.3.7.2	Short run	31
6.3.7.3	Continuous run	32
6.3.8	Softstart and softstop function	32
6.3.9	Rotor selection	33
6.4	Switching the centrifuge off.....	33
7	Malfunctions and error correction.....	34
7.1	General malfunctions.....	34
7.1.1	Emergency lid release.....	35
7.2	Table of error codes	36
7.3	Service contact.....	37

8	Maintenance and service	38
8.1	Maintenance	38
8.1.1	Centrifuge	38
8.1.2	Accessories	39
8.1.2.1	Plastic accessories	39
8.1.3	Rotors	40
8.1.4	Glass breakage	40
8.2	Sterilisation and disinfection of the rotor chamber and accessories	41
8.2.1	Autoclaving	42
8.3	Service	43
8.4	Return of defective units	44
9	Disposal	45
10	Technical data	46
10.1	Ambient conditions	46
11	Appendix	47
11.1	Range of accessories	47
11.1.1	Rotor radii	47
11.2	Speed-gravitational-field-diagram	48
11.3	Table of the service life of rotors and accessories	49
11.4	Resistance data	51
11.5	EC declaration of conformity	55
12	Index	57

1 General information

1.1 Importance of the operating manual

A fundamental requirement for the safe and trouble-free operation of the centrifuge is to be familiar with the fundamental safety instructions and all possible hazards.

The operating manual includes important information concerning the safe operation of the centrifuge.

This operating manual and, in particular, the notes on safety and hazards must be observed by all persons operating the centrifuge.

In addition, the local rules and regulations for the prevention of accidents must be complied with.

1.2 Intended use

Centrifuges are power-driven machines that separate liquids from solid matter, liquid mixtures, or solid mixtures by centrifugal force. They are solely intended for this purpose. Any other use beyond this area of application is regarded as improper use. Sartorius Lab Instruments GmbH & Co. KG cannot be held liable for any damage resulting from such improper use.

The intended use also includes

- observation of all the notes and instructions included in the operating manual and
- compliance with the care, cleaning, and maintenance instructions.

1.3 Warranty and liability

The warranty and liability are subject to our "General Terms and Conditions of Business Relating to Goods and Services of Sartorius" that were distributed to the operator upon the conclusion of the contract.

Warranty and liability claims are excluded if they are due to:

- improper use.
- non-compliance with the safety instructions and hazard warnings in the operating manual.
- improper installation, start-up, operation, or maintenance of the centrifuge.

1.4 Copyright

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

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

Copyright notice:

This operating manual, including all of its components, is protected by copyright. Any use beyond the limits of the copyright law is not permitted without our approval.

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

1.5 Standards and regulations

EC declaration of conformity
(see chapter 11.5 - "EC declaration of conformity")

1.6 Scope of supply

The centrifuge comprises:

- | | |
|--|----------------------------------|
| • 1 connection cable | depending on the voltage variant |
| • 1 spanner, size 5.5
(emergency lid release) | part no. 69CE0001 |
| • 1 rotor fastening nut | part no. 69CE0002 |

Documentation

Operating manual incl. EC declaration of conformity
(see chapter 11.5 - "EC declaration of conformity")

Accessories

according to your order, our order confirmation, and our delivery note.

2 Layout and mode of operation

2.1 Layout of the centrifuge

2.1.1 Functional and operating elements

- 1 Lid
- 2 Display
- 3 User interface
(see chapter 6.3.1 -
"User interface")



Fig. 1: Total view of the centrifuge

- 4 Name plate (see chapter
2.1.2 - "Name plate")
- 5 Mains power switch
- 6 Mains power input



Fig. 2: Rear view of the centrifuge

2.1.2 Name plate

- 1 Manufacturer and registered office
- 2 Type
- 3 Serial number
- 4 Max. density
- 5 Nominal voltage
- 6 Input fuse
- 7 Symbol for special disposal (see chapter 9 - "Disposal")
- 8 CE-mark in compliance with the directive 2006/42/EG
- 9 Part number
- 10 Year of manufacture
- 11 Max. speed
- 12 Max. kinetic energy
- 13 Power consumption



Fig. 1: Example of a name plate

2.2 Mode of operation

2.2.1 Centrifugation principle

Centrifugation is a process for the separation of heterogeneous mixtures of substances (suspensions, emulsions, or gas mixtures) into their components. The mixture of substances, which rotates on a circular path, is subject to centripetal acceleration that is several times greater than the gravitational acceleration.

Centrifuges use the mass inertia inside the rotor chamber for separating the substances. Due to their higher inertia, particles or media with a higher density travel outwards. In doing so, they displace the components with a lower density, which in turn travel towards the centre.

The centripetal acceleration of an object inside a centrifuge, as the effect of centripetal force, depends on the distance between the object and the axis of rotation as well as on the angular velocity. It increases linearly as a function of the distance with regard to the axis of rotation and quadratically as a function of the angular velocity. The bigger the radius in the rotor chamber is and the higher the speed is, the higher the centripetal acceleration is. However, the forces acting on the rotor also increase.

2.2.2 Area of application

Depending on the area of application of the centrifuge and also on the particle size, solids content, and volume throughput of the mixture of substances that is to be centrifuged, there are different types of centrifuges.

The areas of application go from household use as a salad spinner or honey separator up to specialised technical applications in the clinical, biological, or biochemical context:

- For numerous clinical examinations, cellular material must be separated from the liquid to be analysed. The normal separation process can be sped up considerably by using laboratory centrifuges.
- In the metal-working industry, centrifuges are used for separating oil from metal cuttings. Dairies use centrifuges in order to separate cow's milk into cream and low-fat milk.
- Particularly big centrifuges are used in the sugar industry for separating the syrup from the crystalline sugar.
- Ultracentrifuges are predominantly used in biology and biochemistry in order to isolate particles, e.g. viruses. They are specifically designed for high speeds up to 500,000 rpm. The rotor moves in a vacuum in order to avoid air friction.

2.2.2.1 Speed, radius, and relative centrifugal force

The acceleration g , which the samples are subject to, can be increased by increasing the radius in the rotor chamber and by increasing the speed. These three parameters are interdependent and linked with each other via the following formula:

$$\text{Relative centrifugal force RCF} = 11.18 \times 10^{-6} \times r \times n^2$$

r = radius in cm

n = speed in rpm

RCF without any dimension

If two values are entered, the third value is determined by way of the stated formula. If, afterwards, the speed or the radius is changed, the resulting relative centrifugal force will be recalculated automatically by the control unit. If the RCF is changed, the speed will be adapted while the specified radius is maintained.

The speed-gravitational-field-diagram provides an overview of the relationship between speed, radius, and RCF (see chapter 11.2 - "Speed-gravitational-field-diagram").

2.2.2.2 Density

The laboratory centrifuge is suitable for the separation of constituents of different densities in mixtures with a maximum density of 1.2 g/cm³. All information concerning the speed of rotors and accessories refers to liquids with a density corresponding to this specification. If the density is above this value, the maximum permissible speed of the centrifuge must be reduced based on the following formula:

$$n = n_{max} \times \sqrt{(1,2 / Rho)}$$

Rho = density in g/cm³

3 Safety

3.1 Marking of the unit

The following symbols are used for all types of centrifuges manufactured by Sartorius:

	Dangerous voltage		On (Power)
	Hot surface		Off (Power)
	Caution! Risk of bruising		Name plate (see chapter 2.1.2 - "Name plate")
	Protective earth (ground)		CE mark in compliance with the directive 2006/42/EC
	Earth (ground)		Do not dispose as part of domestic waste
	Unplug the mains plug		GS mark (tested safety; only for Germany)
	Arrow indicating the direction of rotation		NRTL mark (only for the USA and Canada)
	Label concerning the FCC rules (only for Canada)		RCM mark (only for Australia)
	Label indicating the emergency release system		China RoHS 2 mark (only for China)

i
NOTE

Safety indications on the centrifuge must be kept readable at all times. If necessary, they must be replaced.

i
NOTE

Not all of the symbols/labels are used for this centrifuge type.

3.2 Explanation of the symbols and notes

In this operating manual, the following names and symbols to indicate hazards are used:



DANGER

This symbol stands for a direct hazard to the life and health of persons.

Non-observance of these symbols causes serious health problems up to life-endangering injuries.



DANGER

This symbol stands for a direct hazard to the life and health of persons due to electrical voltage.

Non-observance of these symbols causes serious health problems up to life-endangering injuries.



WARNING

This symbol stands for a potential hazard to the life and health of persons.

Non-observance of these symbols can cause serious health problems up to life-endangering injuries.



CAUTION

This symbol indicates a potentially hazardous situation

Non-observance of these notes can cause minor injuries or damage to property.



NOTE

This symbol indicates important information.

3.3 Responsibility of the operator

The operator is responsible for authorising only qualified personnel to work on the centrifuge (see chapter 3.4 - "Operating personnel").

The areas of responsibility of the personnel concerning the operation, maintenance, and care of the unit must be clearly defined.

The safety-conscious work of the personnel in compliance with the operating manual and the relevant EC and national health and safety regulations as well as with the accident prevention regulations must be checked at regular intervals (e.g. every month).

Under the international rules for health and safety at work, the operator is obliged to:

- take measures in order to prevent all danger to life or health during work.
- ensure that centrifuges are operated properly and entirely as intended (see chapter 1.2 - "Intended use").
- take protective measures against fire and explosion when working with hazardous substances.
- take measures for the safe opening of centrifuges.

3.4 Operating personnel

Persons operating the unit must

- be familiar with the fundamental regulations concerning workplace safety and accident prevention
- have read and understood this operating manual (and in particular the safety sections and warning notes) and confirmed this with their signature.

3.5 Informal safety instructions

- This operating manual is a part of the product.
- The operating manual must be kept at the location of use of the centrifuge. Ensure that it is accessible at all times.
- The operating manual must be handed over to any subsequent owner or operator of the centrifuge.
- Any changes made must be added to the operating manual.
- In addition to the operating manual, the general and local rules and regulations concerning the prevention of accidents and the protection of the environment must also be supplied.
- Safety and danger indications on the centrifuge must be kept readable at all times. If necessary, they must be replaced.

3.6 Safety instructions

3.6.1 Electrical safety

To reduce the risk of electrical shock, the centrifuge uses a three-wire electrical cord and plug to connect the equipment to earth-ground. To preserve this safety feature:



DANGER

- Ensure that the wall socket is properly wired and grounded.
- Check that the mains voltage agrees with the nominal voltage listed on the name plate.
- Do not place vessels containing liquid on the centrifuge lid or within the safety distance of 30 cm around the centrifuge. Spilled liquids may get into the centrifuge and damage electrical or mechanical components.
- Work on the power supply system must only be performed by certified electricians.
- Inspect the electrical equipment of the unit regularly. Defects such as loose or burnt cables must be eliminated immediately.

3.6.2 Mechanical safety

In order to ensure the safe operation of the centrifuge, observe the following:



WARNING

- Do not open the lid when the rotor is in motion!
- Do not reach into the rotor chamber when the rotor is in motion!
- Do not use the centrifuge if it was installed incorrectly.
- Do not use the centrifuge without panels.
- Do not use the centrifuge if the rotors and inserts show signs of corrosion or other defects.
- Only use the centrifuge with rotors and accessories that have been approved by the manufacturer. In case of doubt, contact the manufacturer (see chapter 7.3 - "Service contact").
- Do not hold your fingers between the lid and the housing when closing the lid. Risk of crushing!
- Defective lid relieving devices could cause the centrifuge lid to fall (contact the service department, if necessary). Risk of crushing!
- Do not hit or move the centrifuge during its operation.
- Do not lean against or rest on the centrifuge during its operation.
- Do not spin any substances that could damage the material of the rotors and buckets of the centrifuge in any way. Highly corrosive substances, for example, damage the material and affect the mechanical strength of the rotors and buckets.
- Stop the centrifuge immediately in the event of a malfunction. Eliminate the malfunction (see chapter 7 - "Malfunctions and error correction") or inform the service department of the manufacturer (see chapter 7.3 - "Service contact").
- Ensure that all repairs are performed only by authorised and specialised personnel.

**WARNING**

- Prior to any start-up, check the centrifuge, rotor, and accessories for signs of damage that can be discerned from the outside. Special attention must be paid to all of the rubber parts (e.g. motor cover, lid seal, and adapters) in terms of visible structural changes. Defective parts must be replaced immediately.
- Open the centrifuge when it is not in use so that moisture can evaporate.

3.6.3 Fire prevention

**DANGER**

- Always use fuses with the same type and rating specified.
- Do not spin explosive or inflammable substances.
- Do not use the centrifuge within hazardous locations.

3.6.4 Chemical and biological safety

If pathogenic, toxic, or radioactive samples are intended to be used in the centrifuge, it is in the responsibility of the user to ensure that all necessary safety regulations, guidelines, precautions, and practices are adhered to accordingly.

**DANGER**

- Infectious, toxic, pathogenic, and radioactive substances may only be used in special, certified containment systems with a bio-seal in order to prevent the material from being released.
- Take suitable precautions for your own safety if there is a risk of toxic, radioactive, or pathogenic contamination
- Materials that chemically react with each other with a high level of energy are prohibited.

**WARNING**

- Keep informed about local measures to avoid harmful emissions (depending on the substances to be centrifuged).
- Protective clothing is not required for the operation of the centrifuge. The materials to be centrifuged may, however, require special safety measures (e.g. centrifugation of infectious, toxic, radioactive, or pathogenic substances).

3.6.5 Safety instructions for centrifugation

For safe operation, observe the following before starting the centrifuge:



WARNING

- Ensure that the centrifuge was set up properly (see chapter 5 - "Set-up and connection").
- Maintain a safety distance of at least 30 cm (12 inches) around the centrifuge.
- Do not store any dangerous goods in the centrifuge area.
- Do not stay in the safety area longer than what is absolutely necessary for the operation of the centrifuge.
- Only use the centrifuge with rotors and accessories that have been approved by the manufacturer. We explicitly warn against the use of equipment of poor quality. Breaking glass or bursting vessels can cause dangerous imbalances at high speeds
- Ensure that rotor and buckets are correctly fitted (see chapter 6.2.2.1 - "Installation of the rotor").
- Observe the instructions on the installation of accessories (see chapter 6.2.2.2 - "Installation of accessories").
- The rotor must be loaded axial symmetrically at equal weights.
- If liquids with a density $> 1.2 \text{ g/cm}^3$ are used, reduce the speed (see chapter 2.2.2.2 - "Density").
- Do not use the centrifuge if the rotor is loaded asymmetrically.
- Do not use the centrifuge with tubes that are excessively long.



WARNING

3.6.6 Resistance of plastics

Chemical influences have a strong effect on the polymeric chains of plastics, and, therefore, on their physical properties. Plastic parts can be damaged if solvents, acids, or alkaline solutions are used.



NOTE

- Refer to the resistance data (see chapter 11.4 - "Resistance data")!

3.6.7 Service life of rotors and accessories

The rotors and accessories have a limited service life.



WARNING

- Perform regular checks (at least once per month) for safety reasons!
- Pay special attention to changes, such as corrosion, cracks, material abrasion, etc.
- After 10 years, they must be inspected by the manufacturer.
- After 50,000 cycles, the rotor must be scrapped for reasons of safety.
- If other data concerning the service life are engraved on the rotor or bucket, these data shall apply accordingly. For example, a bucket with the engraving "max. cycles = 10,000" has a service life of 10,000 cycles, and a rotor with the engraving "Exp. date 02/20" must be scrapped in February 2020 at the latest.



NOTE

- Refer to the table of rotors and accessories with a different service life (see chapter 11.3 - "Table of the service life of rotors and accessories")!

3.7 Safety devices

3.7.1 Lid lock device

The centrifuge can only be started when the lid is properly closed. The electrical lock must be locked. The lid can only be opened when the rotor has stopped. If the lid is opened by way of the emergency release system during operation, the centrifuge will immediately switch off and decelerate brakeless. If the lid is open, the drive is completely separated from the mains power supply, i.e. the centrifuge cannot be started (see chapter 7.1.1 - "Emergency lid release").

3.7.2 System check

An internal system check monitors the data transfer and sensor signals with regard to plausibility. Malfunctions are indicated by error messages with a number in the speed/rcf display (see chapter 7.2 - "Table of error codes").

3.8 Measures in the event of hazards and accidents



DANGER

- If an emergency arises, switch off the centrifuge immediately!
- If in doubt, call the emergency doctor!

3.9 Remaining hazards

The unit was built state- of- the- art and according to the accepted safety rules. However, danger to life and limb of the operator, or of third parties, or impairments of the unit or other material assets cannot be completely excluded when the unit is being used.

- Use the unit only for the purpose that it was originally intended for (see chapter 1.2 - "Intended use").
- Use the unit only if it is in a perfect running state.
- Immediately eliminate any problems that can affect safety.

4 Storage and transport

4.1 Dimensions and weight

	Centrisart® A-14
Height:	176 mm
Height with open lid:	310 mm
Width:	212 mm
Depth:	266 mm
Weight:	6 kg

4.2 Storage conditions

The centrifuge can be stored in its original packaging for up to a year.

- Store the centrifuge only in dry rooms.
- The permissible storage temperature is between -20°C and +60°C.
- If you would like to store it for more than one year, or if you intend to ship it overseas, please contact the manufacturer.

4.3 Notes on transport

- The centrifuge is not equipped with a transport safety device.
- When lifting the centrifuge, always reach under the centrifuge from the side.



CAUTION

The centrifuge weighs approx. 6 kg!

- For transport use suitable packaging and, if at all possible, the original packaging. (see chapter 4.4 - "Packaging").

4.4 Packaging

The centrifuge is packaged in a cardboard box.

- Open the box.
- Take out the box containing the accessories.
- Lift the centrifuge with both foam cushions out of the cardboard box. When lifting the centrifuge, always reach under the centrifuge from the side.



CAUTION

The centrifuge weighs approx. 6 kg!

- Retain the packaging for any possible future transport of the centrifuge.

4.5 Transport safety device

The centrifuge is not equipped with a transport safety device.

5 Set-up and connection

5.1 Installation site

Operate the centrifuge only in closed and dry rooms.

All the energy supplied to the centrifuge is converted into heat and emitted to the ambient air.

- Ensure sufficient ventilation.
- Keep a safety distance of at least 30 cm around the centrifuge so that the vents in the centrifuge remain fully effective.
- Do not subject the centrifuge to thermal stress, e.g. by positioning it near heat generators.
- Avoid direct sunlight (UV radiation).
- The table must be stable and have a solid, even surface.
- Attention: During transport from cold to warmer places, condensational water will collect inside the centrifuge. It is important to allow sufficient time for drying (min. 24 h) before the centrifuge can be used again.

5.2 Power supply

5.2.1 Type of connection



DANGER

The operating voltage on the name plate must correspond to the local supply voltage!

Sartorius centrifuges are units of protection class I. The centrifuges of this model series have a three-wire power cord with an IEC C13 connector.

5.2.2 Customer-provided fuses

Typically, the centrifuge must be protected with 16 Amp L or B fuses that are to be provided by the customer.

6 Using the centrifuge

6.1 Initial start-up



DANGER

- Before the initial start-up, please ensure that your centrifuge is properly set up and installed (see chapter 5 - "Set-up and connection").

6.2 Switching the centrifuge on

- Press the mains power switch.
The display then illuminates. The centrifuge is ready for operation.

6.2.1 Opening and closing the lid

The lid can be opened if the centrifuge is at a standstill.

- Press the lid key in order to open the lid.
The centrifuge cannot be started if the lid is opened.
- To close, press with both hands slightly on the lid until the lid lock is locked.



WARNING

Do not place your fingers between the lid and the housing when closing the lid.
Risk of crushing!

6.2.2 Installation of rotors and accessories

6.2.2.1 Installation of the rotor

- 1 Safety lock
- 2 Snap-on lid
- 3 Rotor fastening nut
- 4 Motor shaft
- 5 Angle rotor

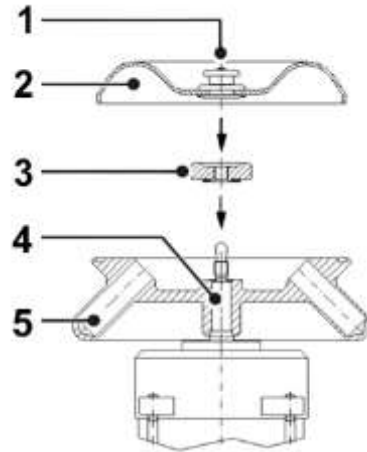


Fig. 2: Installation of rotor and rotor lid

- Open the centrifuge lid by pressing the lid key.
- Unscrew the rotor fastening nut from the motor shaft (counter-clockwise).
- Lower the rotor with its central bore straight down onto the motor shaft.
- Hand-tighten the rotor fastening nut clockwise (corresponds to 1 Nm). The rotor fastening nut and conical spring washer must face the rotor (i.e. the writing must be shown on top).

- 1 Conical spring washer

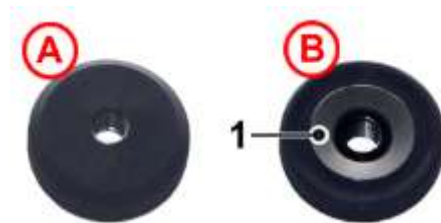


Fig. 3: Rotor fastening nut from above (A) and from below (B)

- To lock the snap-on lid, pull the safety lock in the middle of the lock upwards and put the lid on the rotor. The snap-on lid will be locked when the safety lock is released.



WARNING

Once a day or after 20 cycles, the rotor tie-down screw must be loosened by some turns, and the rotor must be lifted and fastened again. This ensures a proper connection between the rotor and the motor shaft.

- Rotors can be used without a cover. This leads, however, to higher levels of noise and temperatures when running.
- Follow the safety instructions and hazard warnings (see chapter 3 - "Safety")!

6.2.2.2 Installation of accessories

- Only use tubes that are suitable for the rotor (see chapter 11.1 - "Range of accessories").
- Always load the axial symmetrical inserts of the rotors with the same accessories and fill to avoid imbalance.

Centrifugation with low capacity

- Install the tubes axial symmetrically so that the rotor is loaded evenly.
- It is not permissible to load angle rotors on only one axis.

correct



incorrect



Fig. 4: Permissible and impermissible loading of an angle rotor (example illustration)

6.2.2.3 Tubes

- Load the tubes outside of the centrifuge. Liquids in the bores of the rotor cause corrosion.
- Fill the tubes carefully and arrange them according to their weight. Imbalances result in the excessive wear of the bearings.
- Follow the safety instructions and hazard warnings (see chapter 3 - "Safety")!

6.3 Control System "Spincontrol Basic"

6.3.1 User interface

- 1 Display
- 2 Select key
- 3 Arrow keys
- 4 Lid key
- 5 Start/Stop key



Fig. 5: User interface of the centrifuge

The centrifuge is started directly via the user interface. When the centrifuge is switched on, all segments will be illuminated for a short time. It is now ready for operation.

6.3.2 Display

The centrifuge display has the following display fields:

- 1 Speed / RCF field
- 2 Field for acceleration / deceleration curves and run mode
- 3 Time field

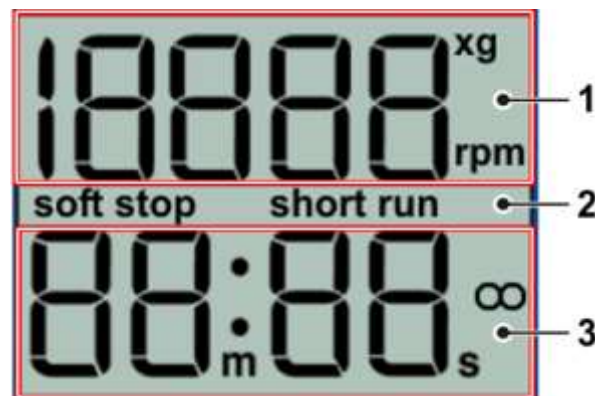


Fig. 6: Display, completely active

6.3.3 Starting a centrifugation run

The centrifuge is ready for operation when the mains power switch is on and the lid is closed.

- Press the start/stop key in order to start a centrifugation run.

6.3.4 Interrupting a centrifugation run

- Press the start/stop key in order to interrupt a centrifugation run. The centrifugation run will be terminated prematurely.

6.3.5 Interrupting a deceleration process

- Press the start/stop key during a deceleration process in order to interrupt it and to restart the centrifuge.

6.3.6 Speed / Relative centrifugal force (RCF)

The RCF value is determined by the rotor geometry and speed. The RCF and speed values, therefore, depend on each other. If one of the two values is entered, the other value will be set automatically.

- To preselect a speed value or a RCF value, press the select key repeatedly until the corresponding unit flashes on the display.
- Select the desired speed or RCF value via the arrow keys.
- Confirm the selection by pressing the start/stop key.

The data will also be accepted and saved after approx. 20 seconds if no other key is pressed within this time period.

During operation, you can switch from the speed value to the RCF value and vice versa via the arrow keys.

6.3.6.1 Changing the speed/RCF value during centrifugation

The preselected speed or RCF value can be changed during centrifugation.

- Press the select key repeatedly until the desired unit flashes on the display.
- Change the speed or RCF value by pressing the arrow keys. The parameters will take effect immediately.

6.3.7 Runtime

The preselected total runtime is displayed in the lower line of the display. During centrifugation, the remaining runtime is displayed. The runtime of the centrifuge can be set at one-second-intervals up to 99 minutes and 59 seconds.

- To select a runtime, press the select key repeatedly until the time unit flashes on the display.
- Select the desired runtime by pressing the arrow keys.
- Confirm the selection by pressing the start/stop key.

The data will also be accepted and saved after approx. 20 seconds if no other key is pressed within this time period.

6.3.7.1 Changing the runtime during centrifugation

The preselected runtime can be changed during centrifugation.

- Press the select key repeatedly until the time unit flashes on the display.
- Change the desired runtime by pressing the arrow keys. The parameters will be accepted immediately.



NOTE

If the centrifugation time is changed during the run, the centrifuge will run for the entire new time and will disregard the previous runtime that has already elapsed.

6.3.7.2 Short run

During the short run, the centrifuge accelerates at maximum power until the maximum speed is reached.

- Keep the quick run key pressed during the short run.

The message "short run" and the duration of the short run is displayed. When the quick run key is released, the centrifuge decelerates at maximum power to a standstill.

After the short run, the lid unlocks automatically and the program that was set beforehand is displayed again.

6.3.7.3 Continuous run

During the continuous run, the runtime of the centrifuge is unlimited and must be stopped manually. The centrifuge accelerates during the continuous run until the set speed is reached.

- To start the continuous run, press the select key until the time unit flashes on the display.
- Press the down-key (arrow key) until the display switches from "00:10" to "--:--".

After 99 min 59 sec, any additional runtime will no longer be displayed, but the centrifugation will continue.

- To stop a continuous run, press the start/stop key. The centrifugation will end.
- Enter a runtime. The centrifugation will end after this time.

6.3.8 Softstart and softstop function

The softstart function is used to extend the acceleration time, whereas the softstop function is used to extend the deceleration time. The current combination is shown on the display.

- To activate the softstart and softstop functions, press the select key repeatedly until "Soft-stop/start" flashes in the upper left part of the display.
- Press the arrow keys until "Soft-stop/start on" is displayed. The softstart and softstop functions are now activated.
- Press the select key until "Soft stop on" is displayed. Only the softstop function will be active and the centrifuge will start at normal speed.
- Press the select key until "Soft off" is displayed. The softstart and softstop functions are now deactivated.

The data will be accepted immediately and saved after approx. 20 seconds.

You can also change the settings as described above during a centrifugation run.

6.3.9 Rotor selection

In the delivery status of the centrifuge, the rotor A2B24x1.5mLRVA14 is preselected. If another rotor is installed the configuration must be changed, so that the allowed maximum speed of the rotor can be reached.

- To select another rotor, press the select key and hold it.
- Press the up-key three times until the rotor selection menu is displayed, then release the select key.
- Choose the number of the installed rotor out of the following list, by pressing the arrow keys:

Code	Rotor
4	YCSR-A2B
6	YCSR-A1C

- Press the select key to confirm the input.

The RCF value will be adjusted automatically to the selected rotor.



NOTE

The rotor selection will be saved under the corresponding program number.

6.4 Switching the centrifuge off

- Open the centrifuge when it is not in use so moisture can evaporate.
- Switch the centrifuge off by pressing the mains power switch.

7 Malfunctions and error correction

7.1 General malfunctions

Malfunctions are indicated by error messages with a number in the speed/RCF display.

In the event of a fatal error (e.g. a defective lid lock), a certain safety time will be counted down on the display. During this time, "ERR" and "SAFE" flash alternately on the display. When the time is up, "OFF" will be displayed.



WARNING

Do not switch the centrifuge off until "OFF" is displayed! This is necessary in order to ensure that the rotor is at a complete standstill.

- Eliminate the source of the problem (see tables below).
- Acknowledge the error messages with the lid key.

Type of error	Possible reason	Correction
No indication on the display	No power in the mains supply	Check fuse in the mains supply
	Power cord is not plugged in	Plug in power cord correctly
	Mains power switch off	Switch mains power switch on
	Lid is not closed correctly	Close the lid
Centrifuge cannot be started: The set speed value is displayed in an unchanged manner	Several	Netz aus/ein. Falls sich der Fehler wiederholt, Service verständigen
	The lid lock is not closed correctly	Open and close lid. If the error occurs again, contact service
Centrifuge decelerates during operation and displays an error from 1 to 18 after powering on	Several	Power off/on. If the error occurs again, contact service
Centrifuge decelerates during operation and displays error 19 after powering on	Several	Quit by pressing the lid key
Temperature value cannot be reached (only for refrigerated centrifuges)	Condenser dirty	Contact service

7.1.1 Emergency lid release

In the event of a power failure, it is possible to manually open the centrifuge lid with a socket spanner (part no. 69CE0001, included in the scope of supply). The opening is located in the base plate in the front quarter on the left-hand side.

- Switch the centrifuge off by actuating the mains power switch and by unplugging the mains power plug.
- Pull the centrifuge a bit over the edge of the table and jack it up (e.g. on two wooden logs). Ensure that the centre of gravity of the centrifuge is still located above the table top. If necessary, the unit must be secured by a second person.
- Insert the spanner for the emergency release into the hole (see the illustration, item 1) and attach it to the shaft of the lid lock motor.
- Turn the shaft carefully clockwise with the aid of the wrench until the lid opens.
- Remove the spanner and pull the centrifuge back on the table.



Fig. :7 Opening for the emergency lid release at the bottom of the centrifuge



WARNING

Do not unlock the lid unless the rotor is at a standstill and unless there is no danger to other persons!

If the lid is opened via the emergency lid release during a centrifuge run, the centrifuge will be switched off immediately and decelerate in a brake-less manner. As a general rule, the lock can be opened more easily with the emergency lid release system if the load on it is relieved by exerting slight pressure on the lid.

7.2 Table of error codes

Error no.	Kind of error	Measures	Note
1-9	System error	<ul style="list-style-type: none"> Allow to slow down Power off/on 	All these errors stop the centrifuge or cause it to decelerate brakeless
10-19	Speedometer error	<ul style="list-style-type: none"> Allow to slow down Power off/on 	
20-29	Motor error	<ul style="list-style-type: none"> Power off Ensure ventilation 	
30-39	EEPROM error	<ul style="list-style-type: none"> Allow to slow down Power off/on 	With error 34, 35, and 36, the centrifuge will stop; with error 37 and 38 only an error message will be given
40-45	Temperature error (only for refrigerated centrifuges)	<ul style="list-style-type: none"> Allow to slow down Power off Allow to cool down Provide better ventilation (only air cooled centrifuges) Provide sufficient water throughput (only water cooled centrifuges) 	
46-49	Imbalance error	<ul style="list-style-type: none"> Allow to slow down Power off Eliminate the imbalance 	
50-59	Lid error	<ul style="list-style-type: none"> Press lid key Close lid Remove foreign matter from the opening of the lid lock device 	With error 50 and 51, the centrifuge will stop
60-69	Process error	<ul style="list-style-type: none"> Allow to slow down Power off/on 	With error 60 message "power failure during run", with error 61, the message "stop after power on"
70-79	Communication error	<ul style="list-style-type: none"> Allow to slow down Power off/on 	
80-89	Parameter error	<ul style="list-style-type: none"> Power off Allow to cool down Provide for better ventilation 	With error 83, error message only
90-99	Other errors	<ul style="list-style-type: none"> Check connections Provide sufficient water throughput (only water cooled centrifuges) 	



NOTE

If it is impossible to eliminate the errors, contact the service!
Please find the corresponding contact data for your region on our website
www.sartorius.com.

7.3 Service contact

In the event of queries, malfunctions, or spare part enquiries:

Please find the corresponding service contact data for your region on our website www.sartorius.com.



NOTE

- If you would like to utilise our service, please state the type of your centrifuge and its serial number.

8 Maintenance and service

The centrifuge, rotor, and accessories are subject to high mechanical stress. Thorough maintenance performed by the user extends the service life and prevents premature failure.



CAUTION

If corrosion or other damage occurs due to improper care, the manufacturer cannot be held liable or subject to any warranty claims.

- Use soap water or other water-soluble, mild cleaning agents with a pH value between 6 and 8 for cleaning the centrifuge and accessories (see also).
- Avoid corrosive and aggressive substances.
- Do not use solvents.
- Do not use agents with abrasive particles.
- Do not expose the centrifuge and rotors to intensive UV radiation or thermal stress (e.g. by heat generators).

8.1 Maintenance

8.1.1 Centrifuge

- Unplug the mains power plug before cleaning.
- Carefully remove all liquids, including water and particularly all the solvents, acids, and alkaline solutions from the rotor chamber using a cloth in order to avoid damage to the motor bearings.
- If the centrifuge has been contaminated with toxic, radioactive, or pathogenic substances, clean the rotor chamber immediately with a suitable decontamination agent (depending on the type of contamination).



WARNING

Take suitable precautions for your own safety if there is a risk of toxic, radioactive, or pathogenic contamination.

- Grease the motor shaft slightly after cleaning (grease for load-bearing bolts part no. 69CE0006).

8.1.2 Accessories



CAUTION

For the care of the accessories, special safety measures must be considered as these are measures that will ensure operational safety at the same time!

- Immediately rinse off the rotor, buckets, or accessories under running water if they have come into contact with any liquids that may cause corrosion. Use a brush for test tubes in order to clean the bores of angle rotors. Turn the rotor upside down and allow it to dry completely.
- Clean the accessories outside the centrifuge once a week or preferably after each use. Adapters should be removed, cleaned and dried.



CAUTION

Do not clean the accessories in a dishwasher!

Cleaning in a dishwasher removes the anodised coating; the result is cracking in areas that are subject to stress.

- If the rotors or accessories have been contaminated with toxic, radioactive, or pathogenic substances, clean them immediately with a suitable decontamination agent (depending on the type of contamination). Take suitable precautions for your own safety if there is a risk of toxic, radioactive, or pathogenic contamination.
- Dry the accessories with a soft cloth or in a drying chamber at approx. 50°C.

8.1.2.1 Plastic accessories

The chemical resistance of plastic decreases with rising temperatures (see chapter 11.4 - "Resistance data").

- If solvents, acids, or alkaline solutions have been used, clean the plastic accessories thoroughly.



WARNING

Plastic accessories must not be greased!

8.1.3 Rotors

The rotors are produced with the highest precision, in order to withstand the permanent high stress from high gravitational fields.

Chemical reactions as well as stress-corrosion (combination of oscillating pressure and chemical reaction) can affect or destroy the metals. Barely detectable cracks on the surface can expand and weaken the material without any visible signs.

- Check the material regularly (at least once a month) for
 - cracks
 - visible damage of the surface
 - pressure marks
 - signs of corrosion
 - other changes.
- Check the bores of the rotors and multiple carriers.
- Replace any damaged components immediately for your own safety.

8.1.4 Glass breakage



CAUTION

In the case of glass breakage, immediately remove all glass particles (e.g. with a vacuum cleaner). Replace the rubber cushions since even thorough cleaning will not remove all glass particles.

Glass particles will damage the surface coating (e.g. anodising) of the buckets, which will then lead to corrosion.

Glass particles in the rubber cushions of the buckets will cause glass breakage again.

Glass particles on the pivot bearing of the load-bearing bolts prevent the buckets and carriers from swinging evenly, which will cause an imbalance.

Glass particles in the rotor chamber will cause metal abrasion due to the strong air circulation. This metal dust will not only pollute the rotor chamber, rotor, and materials to be centrifuged but also damage the surfaces of the accessories, rotors, and rotor chamber.

In order to completely remove the glass particles and metal dust from the rotor chamber:

- Grease the upper third of the rotor chamber with e.g. Vaseline.
- Then, let the rotor rotate for a few minutes at a moderate speed (approx. 2000 rpm). The glass and metal particles will now collect at the greased part.
- Remove the grease with the glass and metal particles with a cloth.
- If necessary, repeat this procedure.

8.2 Sterilisation and disinfection of the rotor chamber and accessories

- Use commercially-available disinfectants such as, for example, Sagrotan[®], Buraton[®], or Terralin[®] (available at chemist's shops or drugstores).
- The centrifuge and the accessories consist of various materials. A possible incompatibility must be considered.
- Before using cleaning or decontamination agents that were not recommended by us, contact the manufacturer to ensure that such a procedure will not damage the centrifuge.
- For autoclaving, consider the continuous heat resistance of the individual materials (see chapter 8.2.1 - "Autoclaving").

Please contact us if you have any queries (see chapter 7.3 - "Service contact").



DANGER

If dangerous materials (e.g. infectious and pathogenic substances) are used, the centrifuge and accessories must be disinfected.

8.2.1 Autoclaving

The service life of the accessories essentially depends on the frequency of autoclaving and use.

- Replace the accessories immediately when the parts show changes in colour or structure or in the occurrence of leaks etc.
- During autoclaving, the caps of the tubes must not be screwed on in order to avoid the deformation of the tubes.



NOTE

It cannot be excluded that plastic parts, e.g. lids or carriers, may deform during autoclaving.

Accessories	Max. temp. (°C)	Min. time (min)	Max. time (min)	Max. cycles
Aluminium buckets	134-138	3	5	-
Aluminium rotors	134-138	3	5	-
Glass tubes	134-138	3	40	-
Polyallomer / polycarbonate rectangular carriers	115-118	30	40	-
Polyallomer / polycarbonate round carriers	115-118	30	40	-
Polyamide buckets	115-118	30	40	10
Polycarbonate / polyallomer lids for angle rotors	115-118	30	40	20
Polycarbonate tubes	115-118	30	40	20
Polyphenylsulfone (PPSU) caps for buckets	134-138	3	5	100
Polypropylene balance adapter for blood-bag systems	115-118	30	40	n.s.
Polypropylene copolymer tubes	115-121	30	40	20
Polypropylene rectangular carriers	115-118	30	40	-
Polypropylene rotors	115-118	30	40	20
Polypropylene round carriers	115-118	30	40	-
Polysulfone caps for buckets	134-138	3	5	100
Polysulfone lids for angle rotors	134-138	3	5	100
Rubber adapters	115-118	30	40	-
Stainless-steel balance weight for blood-bag systems	121	30	30	n.s.
Teflon tubes	134-138	3	5	100

8.3 Service



DANGER

In the event of service work that requires the removal of the panels, there is a risk of electric shock or mechanical injury. Only qualified specialist personnel is authorised to perform this service work.

The centrifuge is subject to high mechanical stress. In order to be able to withstand this high level of stress, high-quality components were used during the production of the centrifuge. Nevertheless, wear cannot be excluded and it may not be visible from the outside. Especially the rubber parts that are – among other things – part of the motor suspension, are subject to ageing.

This is why we recommend having the centrifuge checked by the manufacturer during an inspection once per year in the operating state and once every three years in the dismantled state. Rubber parts should be replaced after three years.

Information and appointments

Please find the corresponding contact data for your region on our website www.sartorius.com.



NOTE

- If you would like to utilise our service, please state the type of your centrifuge and its serial number.

8.4 Return of defective units

Although we exercise great care during the production of our products, it may be necessary to return a unit or accessory to the manufacturer.

In order to ensure the quick and economical processing of returns of centrifuges, spare parts, or accessories, we require complete and extensive information concerning the process. Please fill in the following forms completely, sign them, enclose them with the return package, and send them together with the product.

1. Declaration of decontamination

As a certified company and due to the legal regulations for the protection of our employees and of the environment, we are obliged to certify the harmlessness of all incoming goods. For this purpose, we require a declaration of decontamination.

- The form must be filled in completely and signed by authorised and specialised personnel only.
- Affix the original form in a clearly visible manner to the outside of the packaging.



NOTE

We will return the unit if no declaration of decontamination is provided!

2. Form for the return of defective units

This form is for the product-related data. They facilitate the assignment, and they enable the quick processing of the return. If several units are returned together in one packaging, please enclose a separate problem description for every defective unit.

- A detailed problem description is necessary in order to perform the repair quickly and economically.



NOTE

If the form does not include a description of the malfunction, neither a refund nor a credit note can be issued. In this case, we reserve the right to return the unit to you at your expense.

- Please note on the form if you would like to receive a cost estimate. Cost estimates are only prepared upon request and against charge. If an order is placed, these charges will be offset.

The forms can be downloaded online from www.sartorius.com.

9 Disposal



The packaging is made from environmentally-friendly materials that can be used as secondary raw materials. If you no longer need this packaging, bring it to your local recycling and waste disposal facility according to the regulations applicable in your country. In Germany, you can dispose of this material using the VfW dual system (contract number D-59101-2009-1129). The equipment, including accessories and batteries, must not be disposed of in general household waste, and must be recycled similar to electrical and electronic devices. For further information about disposal and recycling option, please contact your local service staff. The partners listed on the following website can be used for disposals within the EU:

1. Go to <http://www.sartorius.com>.
2. Select the summary under "Service".
3. Then select "Information on Disposal".
4. Adresses for local Sartorius disposal contacts can be found in the pdf-files given on this webpage.



NOTE

Sartorius will not take back equipment contaminated with hazardous materials (ABC contamination) either for repair or disposal.

Service adress for disposal:

Please refer to our website (www.sartorius.com) or contact the Sartorius Service Department for more detailed information regarding repair service adresses or the disposal of your device.

10 Technical data

Manufacturer	Sartorius Lab Instruments GmbH & Co. KG Weender Landstraße 94 - 100 37075 Göttingen Germany
Type:	Centrisart® A-14
<u>Connection requirements</u>	
Electrical connection:	see name plate
Protection class:	I
IP code:	20
Connected load (kVA):	0.157
Power consumption (kW):	0.095
Max. current consumption (A):	0,68 (at 220-240 V / 50-60 Hz) 1,08 (at 100-120 V / 50-60 Hz)
Input fuse (AT):	2.0 (at 220-240 V / 50-60 Hz) 4.0 (at 100-120 V / 50-60 Hz)
<u>Performance data</u>	
Max. speed (rpm):	14,800
Max. capacity (ml):	48
Max. gravitational field (x g)	16,163
Max. kinetic energy (Nm):	1,721
<u>Other parameters</u>	
Time range:	10 sec to 99 min 59 sec, short run, continuous run
<u>Physical data</u>	
Height (mm):	176
Height with opened lid (mm):	310
Width (mm):	212
Depth (mm):	266
Weight (kg):	6.0
Noise level (dB(A)):	<60 (at max. speed)

10.1 Ambient conditions

- The figures are valid for an ambient temperature of $+23^{\circ}\text{C} \pm 2^{\circ}\text{C}$ and a nominal voltage $\pm 10\%$.



NOTE

At a nominal voltage of 100V, a tolerance of +10% / -5% applies.

- For indoor use only.
- Allowable ambient temperature $+5^{\circ}\text{C}$ to $+40^{\circ}\text{C}$.
- Max. relative humidity of air 80% up to 31°C with a linear decrease to 50% relative humidity of air at 40°C .
- Maximum altitude 2,000 m above sea level.

11 Appendix

11.1 Range of accessories

Part-no.	Description	Max. speed (rpm)	Max. gravitational field (x g)
YCSR-A2B	Fixed angle rotor 24 x 1.5/2 ml, for Centrisart® A-14 and A-14C, made of aluminium with plastic lid (polysulfone)	14 800	16 163
YCSR-A1C	Angle rotor for 12 x Vivaspin 500 or 12 x 1.5/2 ml, A-14/A-14C	14 800	15 183

The complete list of accessories can be downloaded from www.sartorius.com.

11.1.1 Rotor radii

The information in the accessories table concerning the radius refers to the values of the respective rotor as shown below. The radius calculation is described in chapter 2.2.2.1 - "Speed, radius, and relative centrifugal force".

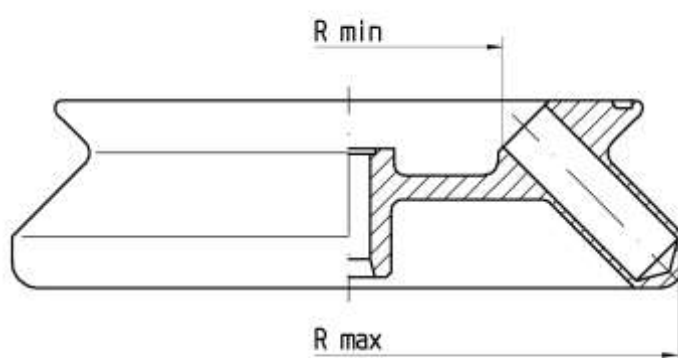


Fig. 8 Minimum and maximum radius of an angle rotor

11.2 Speed-gravitational-field-diagram

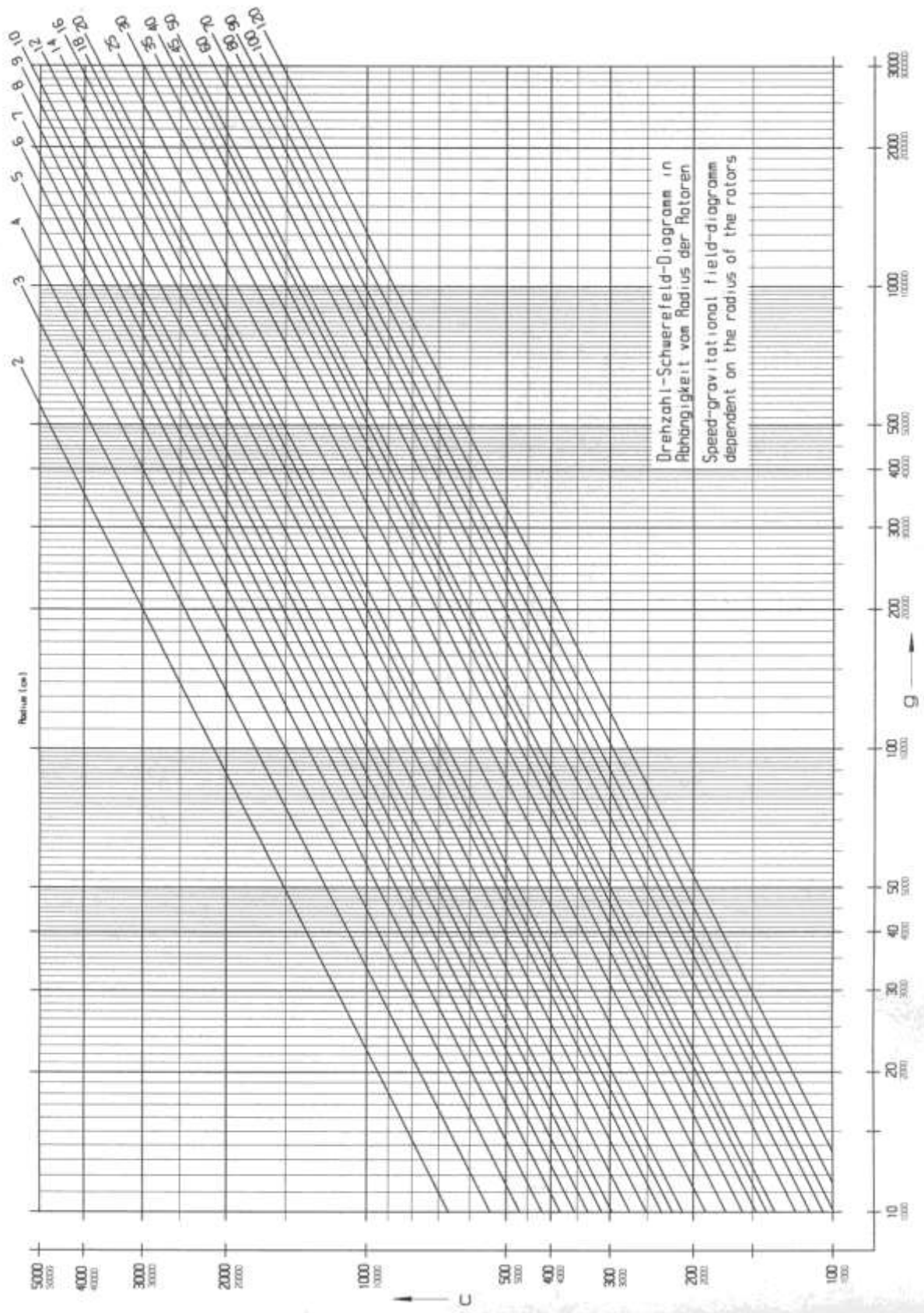



Fig. 9: Speed-gravitational-field-diagram

11.3 Table of the service life of rotors and accessories

Rotors and accessories with a different service life					
If no other data concerning the service life are engraved on the rotor or accessory, rotors and buckets must be checked by the manufacturer after 10 years. After 50,000 cycles, rotors must be scrapped for safety reasons.					
Rotor / Becher	Lebensdauer ("Exp.Date")	Autoklavieren	passend in Zentrifuge		
YCSR-A2B	7 years		Centrisart® A-14, A14C		
YCSR-A1C	5 years	20x	Centrisart® A-14, A14C		

11.4 Resistance data



NOTE

The data refer to resistance at 20°C.

		Concentration	High Density Polyethylene	Polyamide	Polycarbonate	Polyoxymethylene	Polypropylene	Polysulfone	Polyvinyl chloride, hard	Polyvinyl chloride, weak	Polytetrafluorethylene	Acrylonitrile-butadiene-caoutchouc	Aluminium
Medium	Formula	[%]	HDPE	PA	PC	POM	PP	PSU	PVC	PVC	PTFE	NBR	AL
- no data 1 resistant 2 practically resistant 3 partially resistant 4 not resistant													
Acetaldehyde	C ₂ H ₄ O	40	3	2	4	2	3	4	4	-	1	4	1
Acetamide	C ₂ H ₅ NO	saturated	1	1	4	1	1	4	4	-	1	-	1
Acetone	C ₃ H ₆ O	100	1	1	4	1	1	4	4	-	1	4	1
Acrylonitrile	C ₃ H ₃ N	100	1	1	4	3	3	4	4	4	1	4	1
Allyl alcohol	C ₃ H ₆ O	96	1	3	3	2	2	2	2	4	1	1	1
Aluminium chloride	AlCl ₃	saturated	1	3	2	4	1	-	1	-	1	1	4
Aluminium sulfate	Al ₂ (SO ₄) ₃	10	1	1	1	3	1	1	1	1	1	1	1
Ammonium chloride	(NH ₄)Cl	aqueous	1	1	1	2	1	1	1	1	1	1	3
Ammonium hydroxide	NH ₃ + H ₂ O	30	1	3	4	1	1	2	1	-	1	-	1
Aniline	C ₆ H ₇ N	100	1	3	4	1	2	4	4	4	1	4	1
Anisole	C ₇ H ₈ O	100	3	4	4	1	4	4	2	-	1	4	1
Antimony trichloride	SbCl ₃	90	1	4	1	4	1	-	1	-	1	-	4
Benzaldehyde	C ₇ H ₆ O	100	1	3	4	1	1	3	4	4	1	4	1
Benzene	C ₆ H ₆	100	3	2	4	1	3	4	4	-	1	4	1
Boric acid	H ₃ BO ₃	aqueous	1	3	1	2	1	-	-	-	1	1	1
Butyl acrylate	C ₇ H ₁₂ O ₂	100	1	2	4	2	3	4	4	4	1	-	1
Butyl alcohol, normal	C ₄ H ₁₀ O	100	1	1	2	1	1	2	2	4	1	1	1
Calcium chloride	CaCl ₂	alcoholic	1	4	2	3	1	-	-	4	1	1	3
Carbon disulfide	CS ₂	100	4	3	4	2	4	4	4	4	1	3	1
Carbon tetrachloride (TETRA)	CCl ₄	100	4	4	4	2	4	4	4	4	1	3	1
Chlorine	Cl ₂	100	4	4	4	4	4	4	4	4	1	-	3
Chlorine water	Cl ₂ x H ₂ O		3	4	4	4	3	-	3	3	1	-	4
Chlorobenzene	C ₆ H ₅ Cl	100	3	4	4	1	3	4	4	4	1	4	1
Chloroform	CHCl ₃	100	3	3	4	4	3	4	4	4	1	4	3

11 Appendix



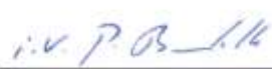
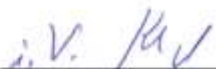
Medium	Formula	Concentration [%]	- no data 1 resistant 2 practically resistant 3 partially resistant 4 not resistant										
			High Density Polyethylene	Polyamide	Polycarbonate	Polyoxymethylene	Polypropylene	Polysulfone	Polyvinyl chloride, hard	Polyvinyl chloride, weak	Polytetrafluorethylene	Acrylonitrile-butadiene-caoutchouc	Aluminium
Medium	Formula	[%]	HDPE	PA	PC	POM	PP	PSU	PVC	PVC	PTFE	NBR	AL
Chromic acid	CrO ₃	10	1	4	2	4	1	4	1	-	1	4	1
Chromic potassium sulphate	KCr(SO ₄) ₂ x 12H ₂ O	saturated	1	2	1	3	1	-	1	-	1	-	3
Citric acid	C ₆ H ₈ O ₇	10	1	1	1	2	1	1	1	1	1	1	1
Citric acid	C ₆ H ₈ O ₇	50	1	3	1	2	1	-	-	-	1	1	1
Copper sulphate	CuSO ₄ x 5H ₂ O	10	1	1	1	1	1	1	1	1	1	1	4
Cyclohexanol	C ₆ H ₁₂ O	100	1	1	3	1	1	1	1	4	1	2	1
Decane	C ₁₀ H ₂₂	100	-	1	2	1	3	-	-	-	1	2	1
Diaminoethane	C ₂ H ₈ N ₂	100	1	1	3	1	1	-	3	4	1	1	1
Diesel fuel	—	100	1	1	3	1	1	-	1	3	1	1	1
Dimethyl formamide (DMF)	C ₃ D ₇ NO	100	1	1	4	1	1	4	3	-	1	3	1
Dimethyl sulfoxide (DMSO)	C ₂ H ₆ SO	100	1	2	4	1	1	4	4	-	1	-	1
Dimethylaniline	C ₈ H ₁₁ N	100	-	3	4	2	4	-	-	-	1	-	1
Dioxane	C ₄ H ₈ O ₂	100	2	1	4	1	3	2	3	4	1	3	1
Dipropylene glycol (mono)methyl ether	C ₄ H ₁₀ O	100	3	1	4	1	4	4	4	4	1	-	1
Ethyl acetate	C ₄ H ₈ O ₂	100	1	1	4	1	1	4	4	4	1	4	1
Ethylene chloride	C ₂ H ₄ Cl ₂	100	3	3	4	1	3	4	4	4	1	-	1
Ferrous chloride	FeCl ₂	saturated	1	3	1	3	1	1	1	1	1	-	4
Formaldehyde solution	CH ₂ O	30	1	3	1	1	1	-	-	-	1	2	1
Formic acid	CH ₂ O ₂	100	1	4	3	4	1	3	3	1	1	2	1
Furfural	C ₅ H ₄ O ₂	100	1	3	3	2	4	-	-	-	1	4	1
Gasoline	C ₅ H ₁₂ - C ₁₂ H ₂₆	100	2	1	3	1	3	3	2	-	1	1	1
Glycerol	C ₃ H ₈ O ₃	100	1	1	3	1	1	1	1	2	1	1	1
Heptane, normal	C ₇ H ₁₆	100	2	1	1	1	2	1	2	4	1	1	1
Hexane, n-	C ₆ H ₁₄	100	2	1	2	1	2	1	2	4	1	1	1
Hydrogen chloride	HCl	5	1	4	1	4	1	1	1	-	1	2	4
Hydrogen chloride	HCl	concentrated	1	4	4	4	1	1	2	3	1	4	4
Hydrogen peroxide	H ₂ O ₂	3	1	3	1	1	1	1	1	-	1	3	3
Hydrogen peroxide	H ₂ O ₂	30	1	4	1	4	1	1	1	-	1	3	3
Hydrogen sulphide	H ₂ S	10	1	1	1	1	1	1	1	3	1	3	1

- no data 1 resistant 2 practically resistant 3 partially resistant 4 not resistant		Concentration	High Density Polyethylene	Polyamide	Polycarbonate	Polyoxymethylene	Polypropylene	Polysulfone	Polyvinyl chloride, hard	Polyvinyl chloride, weak	Polytetrafluorethylene	Acrylonitrile-butadiene-caoutchouc	Aluminium
			Medium	Formula	[%]	HDPE	PA	PC	POM	PP	PSU	PVC	PVC
	Iodine, tincture of	I ₂	1	4	3	1	1	-	4	4	1	1	1
	Isopropyl alcohol	C ₃ H ₈ O	100	1	1	1	1	1	1	4	1	-	2
	Lactic acid	C ₃ H ₆ O ₃	3	1	3	1	2	1	1	2	-	1	1
	Magnesium chloride	MgCl ₂	10	1	1	1	1	1	1	1	1	1	1
	Mercuric chloride	HgCl ₂	10	1	4	1	3	1	1	1	1	1	4
	Mercury	Hg	100	1	1	1	1	1	1	3	1	1	3
	Methyl acetate	C ₃ H ₆ O ₂	100	1	1	4	2	1	-	4	4	1	1
	Methyl alcohol	CH ₄ O	100	1	2	4	1	1	3	1	3	1	2
	Methyl benzene	C ₇ H ₈	100	3	1	4	1	3	4	4	4	1	4
	Methyl ethyl ketone (MEK)	C ₄ H ₈ O	100	1	1	4	1	1	4	4	4	1	4
	Methylene chloride	CH ₂ Cl ₂	100	4	3	4	3	3	4	4	4	1	-
	Mineral oil	—	100	1	1	1	1	1	1	1	-	1	1
	Nitric acid	HNO ₃	10	1	4	1	4	1	1	1	-	1	4
	Nitric acid	HNO ₃	100	4	4	4	4	4	-	4	-	1	4
	Nitrobenzene	C ₆ H ₅ NO ₂	100	3	4	4	3	2	4	4	4	1	4
	Oleic acid	C ₁₈ H ₃₄ O ₂	100	1	1	1	2	1	-	1	-	1	3
	Oxalic acid	C ₂ H ₂ O ₄ x 2H ₂ O	100	1	3	1	4	1	1	1	1	1	2
	Ozone	O ₃	100	3	4	1	4	3	1	1	-	1	4
	Petroleum	—	100	1	1	3	1	1	1	1	3	1	1
	Phenol	C ₆ H ₆ O	10	1	4	4	4	1	4	1	3	1	3
	Phenol	C ₆ H ₆ O	100	2	4	4	4	1	3	4	3	1	3
	Phosphoric acid	H ₃ PO ₄	20	1	4	2	4	1	-	-	-	1	2
	Phosphorus pentachloride	PCl ₅	100	-	4	4	4	1	-	4	4	1	-
	Potassium hydrogen carbonate	CHKO ₃	saturated	1	1	2	1	1	-	-	-	1	-
	Potassium hydroxide	KOH	30	1	1	4	3	1	1	1	1	1	-
	Potassium hydroxide	KOH	50	1	1	4	3	1	1	1	1	1	-
	Potassium nitrate	KNO ₃	10	1	1	1	1	1	-	-	-	1	1
	Potassium permanganate	KMnO ₄	100	1	4	1	1	1	-	1	-	1	3
	Pyridine	C ₅ H ₅ N	100	1	1	4	1	3	4	4	4	1	4
	Resorcinol	C ₆ H ₆ O ₂	5	1	4	2	3	1	4	2	-	1	-
	Silver nitrate	AgNO ₃	100	1	1	1	1	1	1	1	1	1	2

11 Appendix

Medium	Formula	Concentration [%]	- no data 1 resistant 2 practically resistant 3 partially resistant 4 not resistant										
			High Density Polyethylene	Polyamide	Polycarbonate	Polyoxymethylene	Polypropylene	Polysulfone	Polyvinyl chloride, hard	Polyvinyl chloride, weak	Polytetrafluorethylene	Acrylonitrile-butadiene-caoutchouc	Aluminium
Medium	Formula	[%]	HDPE	PA	PC	POM	PP	PSU	PVC	PVC	PTFE	NBR	AL
Sodium bisulphite	NaHSO ₃	10	1	1	2	4	1	-	-	-	1	1	1
Sodium carbonate	Na ₂ CO ₃	10	1	1	1	1	1	-	-	-	1	-	3
Sodium chloride	NaCl	30	1	1	1	1	1	1	1	1	1	1	3
Sodium hydroxide	NaOH	30	1	1	4	1	1	1	1	1	1	2	4
Sodium hydroxide	NaOH	50	1	1	4	1	1	1	1	-	1	2	4
Sodium sulfate	Na ₂ SO ₄	10	1	1	1	1	1	1	1	1	1	1	1
Spirits	C ₂ H ₆ O	96	1	1	1	1	1	1	1	3	1	-	1
Styrene	C ₈ H ₈	100	4	1	4	1	3	-	4	4	1	4	1
Sulphuric acid	H ₂ SO ₄	6	1	4	1	4	1	1	1	-	1	2	3
Sulphuric acid	H ₂ SO ₄	fuming	4	4	4	4	4	4	4	4	1	4	3
Tallow	-	100	1	1	1	1	1	-	1	1	1	1	1
Tetrahydrofuran (THF)	C ₄ H ₈ O	100	3	1	4	1	3	4	4	4	1	3	1
Tetrahydronaphthalene	C ₁₀ H ₁₂	100	3	1	4	1	4	4	4	4	1	-	1
Thionyl chloride	Cl ₂ SO	100	4	4	4	2	4	4	4	4	1	-	3
Tin chloride	SnCl ₂	10	1	4	2	2	1	-	-	-	1	1	4
Transformer oil	-	100	1	1	3	3	1	1	1	-	1	1	1
Trichloroethane	C ₂ H ₃ Cl ₃	100	3	3	4	2	4	4	4	4	1	4	4
Urea	CH ₄ N ₂ O	10	1	1	1	1	1	-	-	-	1	1	1
Urine	-	100	1	1	1	1	1	-	1	1	1	-	2
Vinegar	C ₂ H ₄ O ₂	10	1	4	1	1	1	1	1	1	1	2	1
Vinegar	C ₂ H ₄ O ₂	90	1	4	4	4	1	3	1	4	1	-	1
Wax	-	100	-	1	1		1	-	-	-	1	-	1
Wines	-	100	1	1	1	2	1	1	1	1	1	-	4
Xylene	C ₈ H ₁₀	100	3	1	4	1	4	4	4	4	1	4	1

11.5 EC declaration of conformity

Original		
	EG-/EU-Konformitätserklärung EC / EU Declaration of Conformity	
Hersteller <i>Manufacturer</i>	Sartorius Lab Instruments GmbH & Co. KG 37070 Goettingen, Germany	
	erklärt in alleiniger Verantwortung, dass das Betriebsmittel <i>declares under sole responsibility that the equipment</i>	
Geräteart <i>Device type</i>	Zentrifuge <i>Centrifuge</i>	
Baureihe <i>Type series</i>	Centrisart® A-14	
Modell Model	A-14-1EU, A-14-1US, A-14C-1EU, A-14C-1US	
	in der von uns in Verkehr gebrachten Ausführung allen einschlägigen Bestimmungen der folgenden Europäischen Richtlinien - einschließlich deren zum Zeitpunkt der Erklärung geltenden Änderungen - entspricht und die anwendbaren Anforderungen folgender harmonisierter Europäischer Normen erfüllt: <i>in the form as delivered fulfils all the relevant provisions of the following European Directives - including any amendments valid at the time this declaration was signed - and meets the applicable requirements of the harmonized European Standards listed below:</i>	
2014/30/EU	Elektromagnetische Verträglichkeit <i>Electromagnetic compatibility</i> EN 61326-1:2013	
2011/65/EU	Beschränkung der Verwendung bestimmter gefährlicher Stoffe in Elektro- und Elektronikgeräten (RoHS) <i>Restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS)</i> EN 50581:2012	
2006/42/EG 2006/42/EC	Maschinen <i>Machines</i> EN ISO 12100:2010, EN 61010-1:2010	
	Die Person, die bevollmächtigt ist, die technischen Unterlagen zusammenzustellen: <i>The person authorised to compile the technical file:</i>	
	Sartorius Lab Instruments GmbH & Co. KG International Certification Management 37070 Goettingen, Germany	
	Jahreszahl der CE-Kennzeichenvergabe <i>Year of the CE mark assignment:</i> 16	
	Sartorius Lab Instruments GmbH & Co. KG Goettingen, 2016-04-20	
		
	Dr. Reinhard Baumfalk Vice President R&D	Dr. Dieter Klausgrete Head of International Certification Management
	Diese Erklärung bescheinigt die Übereinstimmung mit den genannten EG- und EU-Richtlinien, ist jedoch keine Zusicherung von Eigenschaften. Bei einer mit uns nicht abgestimmten Änderung des Produktes verliert diese Erklärung ihre Gültigkeit. Die Sicherheitshinweise der zugehörigen Produktdokumentation sind zu beachten. <i>This declaration certifies conformity with the above mentioned EC and EU Directives, but does not guarantee product attributes. Unauthorised product modifications make this declaration invalid. The safety information in the associated product documentation must be observed.</i>	
	Doc: 2033957-01 SU115CE005-01.de,en 1 / 1 PMF: 2033956 OP-113_fo1_2015.10.12	

12 Index

A

Acceleration time	
- see also the softstart and softstop function	32
Accessories, cleaning and care	39
Acid	20, 38, 39
Adapters	39
Alkaline solutions	20, 38, 39
Ambient conditions	46
Ambient temperature	46
Anodised coating	39
Application examples	13
Autoclaving	42

C

Capacity	46
CE mark in compliance with the directive 2006/42/EC	15
Centrifugation of infectious, toxic, radioactive, or pathogenic substances	19
Centrifugation principle	13
Centrifugation with low capacity	28
Centrifuge cannot be started	34
Centrifuge decelerates during operation	34
Centrifuge, cleaning and care	38
Centrifuges	
- definition	9
Changes in colour	42
Changes in structure	42
Changing the runtime during centrifugation	31
Changing the Speed/RCF Value during Centrifugation	30
Chemical and biological safety	19
Chemical resistance of plastic	39
Cleaning agents	41
Cleaning the bores of angle rotors	39
Cleaning the centrifuge	38
Closing the lid	18, 26
Communication error	36
Condensation	25
Connected load	46
Connection cable	10
Contamination	19, 39
Continuous heat resistance	41

Continuous run	32
Copyright	10
Corrosion	18, 21, 28, 38, 39, 40
Cracking	39
Cracks	40
Customer-provided fuses	25

D

Damage of the surface	40
Dangerous goods	20
Dangerous materials	41
Deceleration time	
- see also the softstart and softstop function	32
Declaration of conformity	10, 55
Declaration of decontamination	44
Decontamination agent	39, 41
Deformation of tubes	42
Density	12, 20
Depth	46
Different service life of rotors and accessories	21, 49
Dimensions and weight	23
Direct hazard to the life and health	16
Disinfectants	41
Disinfection of the rotor chamber and accessories	41
Disposal	45
Documentation	10

E

EC declaration of conformity	10, 55
EEPROM error	36
Electrical connection	46
Electrical safety	18
Emergency lid release	35
Error correction	34
Error messages	34
Explanation of the symbols and notes	16
Explosive substances	19

F

Fatal error	34
Fire preventions	19
Form for the return of defective units	44

Functional and operating elements.....	11	Mains voltage	18
G		Maintenance.....	38
General malfunctions	34	Malfunction	34
General Terms and Conditions of Business Relating to Goods and Services of Sartorius.....	9	Malfunctions and error correction.....	34
Glass breakage	40	Manufacturer.....	46
Glass particles.....	40	Marking of the unit	15
Gravitational field.....	46	Max. capacity.....	46
Grease for load bearing bolts.....	38	Max. current consumption	46
H		Max. gravitational field	46
Hazard warnings.....	9, 27, 28	Max. speed.....	46
Height.....	46	Measures in the event of hazards and accidents.....	21
Highly corrosive substances.....	18	Mechanical safety	18
I		Mode of operation	13
Imbalance.....	28	Motor error	36
Imbalance error	36	Motor shaft	27, 38
Importance of the operating manual	9	N	
Important information	16	Name plate.....	12
Infectious substances.....	41	No indication on the display.....	34
Inflammable substances.....	19	No power in the mains supply.....	34
Informal safety instructions	17	Noise level.....	46
Initial start-up.....	26	Nominal voltage	12
Input fuse.....	12	Notes on safety and hazards.....	9
Inspection by the manufacturer.....	43	Notes on transport.....	23
Installation of accessories	28	O	
Installation of rotors and accessories	27	Online download of forms.....	44
Installation of the rotor	27	Opening the lid	26
Installation site.....	25	Operating personnel.....	17
Intended use.....	9	Operating voltage.....	25
Interrupting a centrifugation run.....	30	Operational safety.....	39
Interrupting a deceleration process	30	Overseas shipping.....	23
IP code.....	46	P	
K		Packaging.....	24
Kinetic energy	12	Parameter error.....	36
L		Pathogenic substances.....	19, 38, 41
Layout of the centrifuge.....	11	Performance data.....	46
Leaks	42	Plastic accessories, cleaning and care	39
Lid error	36	Potential hazard to the life and health	16
Lid is not closed correctly.....	34	Potentially hazardous situation	16
Lid lock device.....	21	Power consumption.....	12, 46
Lid lock is not closed correctly.....	34	Power cord is not plugged in.....	34
M		Power supply.....	25
Mains power switch off	34	Pressure marks.....	40
		Prevention of accidents.....	9
		Problem description.....	44

Process error.....	36	Softstart and softstop function.....	32
Protection class.....	25, 46	Solvents.....	20, 38, 39
Q		Spanner for emergency lid release.....	10, 35
Quick run.....	31	Spare fuses.....	10
R		Spare part enquiries.....	37
Radioactive substances.....	19, 38	Speed.....	12, 14, 30, 46
Radius.....	14	Speed-gravitational-field-diagram.....	48
Range of accessories.....	47	Speedometer error.....	36
RCF.....	30	Spincontrol Basic.....	29
Relative centrifugal force (RCF).....	14, 30	Standards and regulations.....	10
Remove glass particles and metal dust from the rotor chamber.....	40	Starting a centrifugation run.....	30
Responsibility of the operator.....	17	Sterilisation of the rotor chamber and accessories.....	41
Return of centrifuges, spare parts, or accessories.....	44	Storage and transport.....	23
Return of defective units.....	44	Storage conditions.....	23
Risk of electrical shock.....	18	Stress-corrosion	
Rotor chamber.....	38	see	
Rotor fastening nut.....	10, 27	corrosion.....	40
Rotor radii.....	47	Structural changes.....	19
Rotor selection.....	33	Supply voltage.....	25
Rotors and accessories with a different service life.....	21, 49	Switching the centrifuge off.....	33
Rotors, cleaning and care.....	40	System check.....	21
Runtime.....	31	System error.....	36
S		T	
Safety area.....	20	Table of error codes.....	36
Safety devices.....	21	Table of rotors and accessories with a different service life.....	21, 49
Safety distance.....	18, 20, 25	Technical data.....	46
Safety instructions.....	9, 27, 28	Temperature error.....	36
Safety instructions for centrifugation.....	20	Temperature value cannot be reached.....	34
Safety, chemical and biological.....	19	Thermal stress.....	38
Safety, electrical.....	18	Time range.....	46
Safety, mechanical.....	18	Toxic substances.....	19, 38
Safety-conscious work.....	17	Transport safety device.....	24
Scope of supply.....	10	Type.....	12
Serial number.....	12, 37, 43	Type of connection.....	25
Service.....	43	Type of the centrifuge.....	37, 43, 46
Service contact.....	37	U	
Service life.....	38	Units of protection class I.....	25
Service life of rotors and accessories.....	21, 49	User interface.....	29
Service life of the accessories.....	42	UV radiation.....	25, 38
Service work.....	43	V	
Set-up and connection.....	25	Vents.....	25
Short run.....	31		

W		Width	46
Warranty and liability.....	9	Work on the power supply system	18
Wear	43	Y	
Weight.....	46	Year of manufacture	12

Sartorius Lab Instruments GmbH & Co. KG
Weender Landstrasse 94–108
37075 Goettingen, Germany

Phone: +49.551.308.0
Fax: +49.551.308.3289
www.sartorius.com

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

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

Copyright notice:

This operating manual, including all of its components, is protected by copyright. Any use beyond the limits of the copyright law is not permitted without our approval. This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:
07/2015