



sartorius

## Centrisart® G-26C

Universal High-Speed Centrifuge  
for laboratory applications

### Advantages

- High speed 62,000 × g: Particularly beneficial to separate even smallest particles
- High clarity: Process library with program preview
- Security: Optional multistep program protection
- Volume | Capacity: 6 × 85 ml | 510 ml



### Product Description

The refrigerated centrifuge Centrisart® G-26C stands out through its high speed with up to 26,000 rpm and a gravitational field with up to 62,000 × g. Therefore the Centrisart® G-26C centrifuge is particularly suitable for the sedimentation of very small particles with low density differences and for specific applications like the isolation of viruses, organelles or membranes, and the precipitation of proteins.

Also, the cooling centrifuge Centrisart® G-26C is a perfect solution for applications that require lower rotation speed, such as isolation of nucleic acids using spin-pillars, or ultrafiltration using Vivaspin® units for sample concentration of e.g. antibody solutions, for sample desalting or buffer exchange, and removal of contamination with small or very large molecular weights. Offering a comprehensive selection of rotors and adapters for different vials as well as the variably adjustable speed range between 100 rpm and 26,000 rpm the Centrisart® G-26C provides an extremely high flexibility.

Quick and easy programming of the refrigerated centrifuge Centrisart® G-26C is facilitated by the use of one button only to select and set all relevant parameters. The centrifuge features a process library with tabular program preview and a storage capacity for 60 programs. For convenient management all stored programs are displayed with name and values for rotation speed, run time, and temperature. Searching for, selection of, and even starting a specific program directly from the process library is

exceptionally simple and highly convenient. Optionally, functions like editing, saving, loading, and starting of individual programs can be password protected to prevent unauthorized access. Applying the function "program rotation" several programs can be combined and executed in succession.

The innovative curve editor of the refrigerated high-speed micro-centrifuge Centrisart® G-26C provides the option to individually configure acceleration and deceleration curves in accordance with the specific sample material. Additional protection of the samples is guaranteed through the high-performance compressor which reliably cools the rotor chamber to –20° C. The temperature monitoring function only tolerates runs to be started when the temperature is within the permitted tolerance range. During the run the cooling system ensures reliable sample cooling to 4°C even at maximum rotation speed. The integrated run monitoring function continuously controls the rotation speed and run time during a centrifugation process, compares the actual values to the programmed values and displays a compact summary after each run.

## Additional Advantages

- Optimum combination: Rotors for Sartorius VivaSpin® filtration units
- Smart Control: Setting all run parameters using just one control knob
- Sample protection: Standstill cooling and “Rapid temp” pre-cooling

## Special Features

- Setting run parameters using just one control knob
- Splash-proof keypad with clear display
- Backlit keys with tactile feedback
- Operation with gloves possible
- Automatic conversion of speed and RCF
- Width only 57 cm
- Speed 100 – 26,000 rpm, adjustable in steps of 100 rpm
- Max. gravitational field (RCF) 61,973 g, adjustable in increments of 10 × g
- Time 10 sec. – 99 min., adjustable in steps of 1 sec., short run, continuous run
- Optional runtime counting after reaching the set speed
- Continuous run monitoring
- Automatic rotor and imbalance detection
- Cycle counter
- Optional soft start or soft stop adjustable in 10 steps, configurable
- Process library for 60 programs
- Optional access protection for program functions
- Reliable sample cooling to 4°C even at max. rpm
- “Rapid temp” precooling
- Stand still cooling
- Maintenance-free motor
- Two motorized lid locks, easy lid opening supported by two pneumatic springs
- Emergency lid lock release with no need to open the casing to access
- High chemical resistance
- Pleasantly quiet operation

## Technical Specifications

Max. speed rpm	26,000
Max. capacity	6 × 85 ml
Max. gravitational field	61,973 g
Min. speed rpm	100
Temperature range	– 20°C to + 40°C
Temperature increments	Adjustable in increments of 1°C
Quick cooling	Quick cool program, rotor cooling even possible during downtimes
Coolant	CFC-free (R404a)
Display	LCD, 116 × 87 m, backlit keys
Control panel	Easy, intuitive 1-button operation
Motor	Maintenance-free, brushless drive motor
Signal	Acoustic signal, can be switched off, configurable signal length
Lid-lock release	Motorized locked lid, automatic lid-lock release
Lid	Inspection window for independent speed measurement, easy-to-open lid port
Rotor	Rotor detection for fixed angle rotors, swinging rotors
Rotor chamber	Stainless-steel
Imbalance safety mechanism	Automated imbalance control
Timer	10 sec. – 99 min. 59 sec., short run   continuous run

Programs	60
Acceleration curves	10, linear, quadratic, user-definable
Deceleration curves	10, linear, quadratic, user-definable
Rotor safety	Automatic detection, cycle counting of the rotors
Dimensions (W × D × H)	570 × 616 × 460 mm
Height with open lid	890 mm
Weight	100 kg
Noise level at max. speed	<63 dB
Power consumption	1,800 W

## Order Information

Order Number	Description
G-26C-1EU	Centrisart® G-26C high-speed refrigerated centrifuge, 220–240 V, 50 Hz
G-26C-1US-2P	Centrisart® G-26C high-speed refrigerated centrifuge, 208 V, 60 Hz, 2 phases
G-26C-1US-3P	Centrisart® G-26C high-speed refrigerated centrifuge, 220–240 V, 60 Hz, 3 phases
G-26C-1JP-2P	Centrisart® G-26C high-speed refrigerated centrifuge, 200 V, 50   60 Hz, 2 phases

## Accessories

Order Number	Description
YCSR-A5B	Fixed angle rotor for 6 × 50 ml round bottom culture tubes (Ø 28.8 × 107 mm), incl. hermetically sealing aluminium lid
YCSR-A6B	Fixed angle rotor for 24 × 1.5   2.0 ml conical reaction tubes, incl. hermetically sealing aluminium lid
YCSA-1	Adapter for 0.25   0.4 ml reaction tubes for rotors YCSR-A2B, YCSR-A4B and YCSR-A6B, 1 set = 2 pcs.
YCSA-2	Adapter for 0.5   0.75 ml reaction tubes Ø 7.9/10 × 28/31 mm for rotors YCSR-A2B, YCSR-A4B and YCSR-A6B, 1 set = 2 pcs.
YCSA-4	Adapter for 0.2 ml PCR reaction tubes for rotors YCSR-A2B, YCSR-A4B and YCSR-A6B, 1 set = 2 pcs.
YCSA-3	Adapter for > 20.000 g for 1.5 ml reaction tubes in rotor YCSR-A6B, 1 set = 2 pcs.
YCSR-A7B	Fixed angle rotor for 4 × 80 85 ml round bottom culture tubes (Ø 38 × 112 mm), incl. hermetically sealing aluminium lid
YCSR-A8B	Fixed angle rotor for 6 × 85 ml round bottom culture tubes (Ø 38 × 112 mm), incl. hermetically sealing aluminium lid
YCSR-A0B	Fixed angle rotor for 6 × 50 ml conical culture tubes, incl. hermetically sealing aluminium lid
YCSR-S3B	Swing-out rotor, 4 place, for bucket YCSB-B9B
YCSB-B9B	Round bucket for adapters YCSA-R0B – YCSA-RDB, sealable with lid YCSM-B9B, 1 set = 2 pcs.
YCSA-R9B	Adapter for YCSB-B9B for 1 conical culture tube 50 ml, 1 set = 2 pcs.
YCSA-R0B	Adapter for YCSB-B9B for 2 conical culture tubes 15 ml, lid max. Ø 23 mm, 1 set = 2 pcs.
YCSA-RAB	Adapter for YCSB-B9B for 4 conical reaction tubes 1.5   2.0 ml, max. Ø 11 mm, 1 set = 2 pcs.
YCSA-RBB	Adapter for YCSB-B9B for 1 flat bottom culture tube 30 ml, Ø 25/31 × 65–95 mm, 1 set = 2 pcs.
YCSA-RCB	Adapter for YCSB-B9B for 3 conical reaction tubes 5.0 ml, 1 set = 2 pcs.
YCSA-RDB	Rubber cushion for YCSB-B9B for 100 ml glas tube, 1 set = 2 pcs.
YCSM-B9B	Round polysulfone sealing cap for round bucket YCSB-B9B, clear, 1 set = 2 pcs.
YCST-G26C	Table for centrifuge made of sheet-steel, varnished, with 2 doors and space for accessories, movable on lockable castors w 570, d 616, h 481 mm

## Technical Data Rotors

Order Number	Rmax [mm]	Rmin [mm]	rpm	RCF
YCSR-S3B	142	38	5,000	3,969
YCSR-A5B	84	21	21,000	41,415
YCSR-A6B	82	50	26,000	61,973
YCSR-A7B	91	14	20,000	40,695
YCSR-A8B	98	28	15,500	26,323
YCSR-A0B	93	31	14,000	20,379

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

Phone +49.551.308.0  
Fax +49.551.308.3289  
www.sartorius.com

USA Toll-free +1.800.635.2906  
UK +44.1372.737159  
France +33.1.70.62.50.00  
Italy +39.039.4659.1  
Spain +34.913.586.095  
Russian Federation +7.812.327.53.27  
Japan +81.3.3740.5408

Specifications subject to change without notice.  
Copyright Sartorius Lab Instruments GmbH & Co. KG.  
Printed in the EU on paper bleached without chlorine.  
Publication No.: W--2108-e161106  
Order No.: 98649-019-52  
Ver. 11 | 2016