



## Centrisart® A-14C

### Micro-Centrifuge

#### Advantages

- Optimum combination – Rotors for Sartorius VivaSpin® filtration units
- Sample protection – Reliable sample cooling to 4 °C at max. speed
- Compact – Very small footprint (27 × 24 cm)
- Volume | Capacity: 24 × 2 ml | 48 ml



#### Product Description

The microcentrifuge Centrisart® A-14C is extremely compact, space-saving. The powerful centrifuge provides a maximum speed of 15,000 rpm and a capacity of 24 × 2 ml. Microcentrifuge A-14C is ideal suited for nucleic acid isolation and purification using spin-columns and concentrating biological samples using Sartorius Vivaspin® 500 units.

Processes like sample concentration of e.g. antibody solutions by ultrafiltration, desalting or buffer exchange, as well as removal of contaminations with small or very large molecular weights, are particularly supported through the special Vivaspin® 500 ultrafiltration unit-tailored rotor YCSR-A1C. In addition, the microcentrifuge is well-suited to sediment cells or particles in small volumes. The device provides large control buttons and easy-to-read displays to facilitate quick and easy input or selection of centrifugation parameters and convenient operation.

Thanks to its small size the microcentrifuge Centrisart® A-14C fits on any laboratory bench. The low height and the motorized lid lock ensure convenient sample loading and removal. Upon a gentle push the lid closes automatically. On completion the lid can be unlocked by a single key stroke and lifts autonomously. With just one finger the lid can then be opened completely and is securely kept open by a spring mechanism. During centrifugation runs the electronic lid lock prevents accidental opening of the lid to provide extensive user protection.

Microcentrifuge Centrisart® A-14C represents exceptional product quality and meets the highest safety requirements. Amongst others, the high quality is reflected by the low-level noise of max. 53 dB and the chemical resistance of the materials processed. Spreading of aerosols is safely prevented by a rotor cover with quick-locking function and both, the rotor chamber and the rotors can easily be cleaned. For temperature-sensitive samples, the active cooling option of the microcentrifuge Centrisart® A-14C can be used. The pre-cool function aid the rapid cool down to the set target temperature and even at maximum speed the model A-14C ensures reliable sample cooling to 4 °C.

## Special Features

- Splash-proof keypad with clear display
- Keys with tactile feedback
- Operation with gloves possible
- Automatic conversion of speed and RCF
- Speed 200 – 15,000 rpm, adjustable in steps of 100 rpm
- Max. gravitational field (RCF) 16,602 g, adjustable in increments of 10 x g
- Timer 10 sec. – 99 min., adjustable in steps of 1 sec., short run, continuous run
- Optional soft start or soft stop
- Memory of the last setting
- Maintenance-free motor
- Motorized lid lock, automatic lid release
- Emergency lid lock release
- Ergonomic operation of the lid with one finger
- High chemical resistance
- Pleasantly quiet operation
- Memory for 10 programs
- Temperature range -10 – + 40°C, adjustable in steps of 1°C
- Reliable sample cooling to 4°C even at max. rpm
- Rapid pre-cooling mode

## Technical Specifications

Max. speed rpm	15,000
Max. capacity	24 × 2 ml
Max. gravitational field	16,602 g
Min. speed rpm	200
Display	LCD, 65 × 47 mm
Motor	Maintenance-free, brushless drive motor
Timer	10 sec. – 99 min. 59 sec., short run   continuous run
Lid-lock release	Motorized locked lid, automatic lid-lock release
Lid	Inspection window for independent speed measurement
Rotor	Fixed angle rotor
Programs	10
Acceleration curves	2
Deceleration curves	2
Dimensions (W × D × H)	273 × 515 × 236 mm
Height with open lid	404 mm
Weight	19 kg
Noise level at max. speed	53 dB
Power consumption	230 W

## Order Information

Order Number	Description
A-14C-1EU	Centrisart® A-14C centrifuge 220 – 240 V; 50   60 Hz
A-14C-1US	Centrisart® A-14C centrifuge 110 – 120 V; 50   60 Hz
A-14C-1EU-SET	Set Centrisart® A-14C centrifuge, 220 – 240 V, 50   60 Hz, incl. rotor YCSR-A1C
A-14C-1US-SET	Set Centrisart® A-14C centrifuge, 100 – 120 V, 50   60 Hz, incl. rotor YCSR-A1C

### Accessories

Order Number	Description
YCSR-A2B	Fixed angle rotor for 24 × 1.5   2.0 ml conical reaction tubes, incl. polysulfone lid YCSM-A1B (not available for A-14-1US)
YCSR-A1C	Fixed angle rotor made of polypropylene for 12 × 1.5   2.0 ml conical reaction tubes or 12 × Vivaspin® 500, incl. polysulfone lid YCSR-A1C
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.
YSCA-4	Adapter for 0.2 ml PCR reaction tubes for rotors YCSR-A2B, YCSR-A4B and YCSR-A6B, 1 set = 2 pcs.

### Technical Data Rotors

Order Number	Rmax [mm]	Rmin [mm]	rpm	RCF
YCSR-A2B	66	30	15,000	16,602
YCSR-A1C	62	29	14,800	15,596

# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

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

Phone +49.551.308.0  
Fax +49.551.308.3289

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan

Phone +33.1.70.62.50.00  
Fax +33.1.64.59.76.39

### Austria

Sartorius Austria GmbH  
Modecenterstrasse 22  
1030 Vienna

Phone +43.1.7965760.0  
Fax +43.1.7965760.24

### Belgium

Sartorius Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles

Phone +32.2.756.06.90  
Fax +32.2.481.84.11

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki

Phone +358.9.755.951  
Fax +358.9.755.95.200

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi

Phone +3623.457.227  
Fax +3623.457.147

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11

Phone +353.1.8089050  
Fax +353.1.8089388

### Italy

Sartorius Italy S.r.l.  
Via Torino 3/5  
20814 Varedo (MB)

Phone +39.039.4659.1  
Fax +39.039.4659.88

### Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001  
Fax +31.30.60.52.917

[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn

Phone +48.61.6473830  
Fax +48.61.6473839

### Russian Federation

LLC "Sartorius RUS"  
Uralskaya str. 4, Lit. B  
199155 St. Petersburg

Phone +7.812.327.53.27  
Fax +7.812.327.53.23

### Spain & Portugal

Sartorius Spain, S.A.  
Avda. de la Industria, 32  
Edificio PAYMA  
28108 Alcobendas (Madrid)

Phone Spain +34.913.586.095  
Phone Portugal +351.800.855.800  
Fax Spain +34.913.589.623  
Fax Portugal +351.800.855.799

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00  
Fax +41.44.746.50.50

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ

Phone +44.1372.737159  
Fax +44.1372.726171

### Ukraine

LLS "Sartorius RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine

Phone +380.44.411.4918  
Fax +380.50.623.3162

## Americas

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
Fax +1.631.254.4253

### Argentina

Sartorius Argentina S.A.  
Int. A. AVALOS 4251  
B1605ECS Munro  
Buenos Aires

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900  
Fax +55.11.4362.8901

### Canada

Sartorius Canada Inc.  
2179 Dunwin Drive #4  
Mississauga, ON L5L 1X2

Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepetzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepetzotlan,  
Estado de México,  
C.P. 54605

Phone +52.55.5562.1102  
Fax +52.55.5562.2942

[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

### Peru

Sartorius Peru S.A.C.  
Av. Emilio Cavenecia 264 San Isidro  
15073 Lima, Perú

Phone +51.1.441 0158  
Fax +51.1.422 6100

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175

Phone +61.3.8762.1800  
Fax +61.3.8762.1828

### China

Sartorius (Shanghai) Trading Co., Ltd.  
3rd Floor, North Wing, Tower 1  
No. 4560 Jinke Road  
Zhangjiang Hi-Tech Park  
Pudong District  
Shanghai 201210, P.R. China

Phone +86.21.6878.2300  
Fax +86.21.6878.2882

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwun Tong  
Kowloon, Hong Kong

Phone +852.2774.2678  
Fax +852.2766.3526

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India

Phone +91.80.4350.5250  
Fax +91.80.4350.5253

### Japan

Sartorius Japan K.K.  
4th Fl., Daiwa Shinagawa North Bldg.  
8-11, Kita-Shinagawa 1-chome  
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408  
Fax +81.3.3740.5406

### Malaysia

Sartorius Malaysia Sdn. Bhd  
Lot L3-E-3B, Enterprise 4  
Technology Park Malaysia  
Bukit Jalil  
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622  
Fax +60.3.8996.0755

### Singapore

Sartorius Singapore Pte. Ltd  
1 Science Park Road,  
The Capricorn, #05-08A,  
Singapore Science Park II  
Singapore 117528

Phone +65.6872.3966  
Fax +65.6778.2494

### South Korea

Sartorius Korea Ltd.  
8th Floor, Solid Space B/D,  
PanGyoYeok-Ro 220, Bundang-Gu  
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700  
Fax +82.31.622.5799

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310

Phone +66.2643.8361-6  
Fax +66.2643.8367



◀ [www.sartorius.com](http://www.sartorius.com)