



BIOSTAT® B-DCU II

O₂ Enrichment



The BIOSTAT® B-DCU II is the second generation of a well proven Fermentor | Bioreactor system, designed for meeting demanding requirements in both research and process development. Unrivalled for scale-down and scale-up modeling of various culture processes, the BIOSTAT® B-DCU II provides a new level of power and flexibility. The BIOSTAT® B-DCU II with integrated O₂-Enrichment gassing capability enables high oxygen transfer for high cell density cultures as well as for sheer-stress sensitive gassing for filamentous organisms. Furthermore, it may help to solve foaming problems due to reduced gassing rates.

Control Tower

- Intuitive touch screen operation
- Industrial controller hardware
- Controller for Agitation, Temperature, pH, pO₂, Foam
- Turbidity, Redox or culture vessel pressure measurement and control
- Automatic gas composition controlled by pO₂-controller
- User configurable 5-stage pO₂ cascade control via agitation speed, O₂-Enrichment with optional MFC's and substrate feeding
- Up to 4 feed controller per vessel
- Controller status indication
- Gas and pump totalizers
- Automatic single and group sensor calibration
- In-process pH-recalibration
- Trend display for up to 8 process values
- Up to 4 balance connections

Supply Tower

"O₂-Enrichment" Gassing System

- Gas mixing of Air and O₂
- O₂-Enrichment capability controlled via DO controller
- Optional mass flow controller for Total Sparger flow or for Air and O₂

Pumps

- 3 integrated pumps, expandable up to 6
- Up to 4 feeding pumps

Temperature System

- Powerful heater (1 kW)
- Integrated controlled cooling water valve
- Circulation pump
- Temperature range 8°C above cooling water up to 80°C

Agitation System

- Speed range 20 up to 2,000 rpm
- Maintenance free
- High torque for power full mixing
- Gear-free for quiet operation

Culture vessel

- Jacketed culture vessel fully equipped with:
- Sensors for Temperature, DO, pH and Foam
 - Stirrer shaft with single mechanical seal
 - Rushton impeller
 - Baffle assembly (not for 0.5 L vessel)
 - Aeration tube with ring Sparger, sterile filters and exhaust cooler
 - Manual sampler with sampling pipe
 - Removable addition bottle support (not for 0.5 L vessel)
 - Addition bottles with stainless steel head piece and sterile filters
 - Inoculation | addition septum port
 - Multiple way addition port
 - Tube, O-ring and tool kit

SCADA Software MFCS/DA

- To accelerate your research activities, a powerful supervisory software MFCS/DA for extended visualization, data acquisition and trend display is included.
- Plug and Play configuration
 - Batch-oriented software package
 - Online data acquisition
 - Sample Data Management
 - Enhanced Plotting
 - Export functions
 - Easy to use programming interface

Features

- Powerful industry rated DCU-4 control system with 15" TFT color touchscreen
- Independent process control for up to six culture vessels
- UniVessel® from 0.5 L to 10 L working volume
- Up to six integrated peristaltic pumps with choices for fixed and analogue speed pumps
- Choice of polarographic or optical pO₂-sensors
- Superior gas mixing with up to 2 Rotameter and Mass Flow Controller
- Culture vessel pressure control
- Easy on-site Supply Tower | Culture vessel upgrade
- Improved connectivity of utilities and probes
- Inclusive Supervisory Process control software
- Validation support available, inclusive Logbook and 3-Level password protection

Technical Specification

Space requirements | Environmental conditions

Space requirement 1-fold 2-fold 3-fold 4-fold 5-fold 6-fold [W×H×D] (without options)	800 1200 1700 2050 2550 3000×780 (10L: 820)×800 [mm]
Space requirement autoclave Ø H (with BB-8844593 flexible adaptor for exhaust cooler) 0.5 L 1 L 2 L 5 L 10 L without tray for storage bottles	170×340 (N/A) 240×500 (340) 270×550 (400) 300×700 (510) 350×820 (620) [mm]
Ambient temperature relative humidity (non-condensating)	5–40°C 85%

Control Tower

Housing material	Stainless steel AISI 304
Display	Touch Screen 15"
Resolution	1024×768 dpi
Communication Control Tower/Supply Tower Control Tower/Host	Ethernet Ethernet

Measurement ranges | resolution

Stirrer speed 0.5 L 1 L 2 L 5 L 10 L (200W) 10 L (400 W)	tbd 20–2000 20–2000 20–1500 20–800 20–1200 [rpm] 1 rpm
Temperature	0–150°C 0.1°C
pH	2–12 0.01 pH
pO ₂	0–250% 1%
Foam and Level	on/off 4 user selectable sensitivities
Turbidity (option)	0–6 AU 0.01 AU
Redox (option)	–2000–2000 mV 1 mV
Pressure (option)	0–1000 mbarg 1 mbar

Gassing System "O₂-Enrichment"

Outlet design	Air aeration with O ₂ supplementation
---------------	--

Flowmeter

Gas flow range "Sparger" 0.5 L 1 L 2 L 5 L 10 L	Air calibrated @ 1.21 bara 20 °C
Accuracy	0.11–1.05 0.16–1.6 0.42–4.2 1.3–13 2–20 [l/min]

Thermal Mass Flow Controller (option)

Flow range "Sparger" Total Flow 1 L 2 L 5 L 10 L	Air calibrated
Accuracy	0.02–1.0 0.06–3 0.06–3 0.2–10 0.4–20 [slpm]

Agitation

Motor performance torque 0.5 L 1 L, 2 L, 5 L, 10 L 10 L	Maintenance and gear-free Servo drive
Maximum impeller tip speed 0.5 L 1 L 2 L 5 L 10 L (200 W) 10 L (400 W)	75/0.28 200/1.03 400/1.32 option [W]/[Nm]

Integrated pumps

Pump head	on/off controlled, pulse-width modulated controlled
Rotation speed	Watson Marlow 102R, for tubing with 1.6 mm wall thickness
Flow rate tube dependent (bore×wall tubing) 1.6×1.6 3.2×1.6 [mm]	20 rpm

Integrated feed pump (option)

Pump head	Speed controlled
Rotation speed	Watson Marlow 102R, for tubing with 1.6 mm wall thickness
Flow rate tube dependent (bore×wall tubing) 1.6×1.6 3.2×1.6 [mm]	1–10 5–50 [rpm]

Temperature control system

Temperature control range	Thermostat system with recirculation pump and solenoid valve for cooling water
Electrical heater	8°C above cooling water to 80°C
Connections to culture vessel Exhaust cooler	1000 W

External connections

Balance standard option/interface/connector	Quick couplings Quick couplings
CO ₂ exhaust analyzer interface/connector	2 4/RS232/M12
Feed pump connection qty./interface/connector	1/RS232/M12
External inputs qty./interface/connector	2/0–10 V/M12

Culture vessel

	0.5 L	1 L	2 L	5 L	10 L
Design	Jacketed glass vessel with stainless steel head plate				
Total volume	0.75	1.6	3	6.6	13 [L]
Working volume	0.15–0.5L	0.4–1	0.6–2	0.6–5	1.5–10 [L]
Headplate ports 19 mm 12 mm 6 mm	– 6 4	3 2 6	3 2 9	3 3 8	7 2 9
Volume storage bottles	250	250	250	500	500 [mL]
pO ₂ sensor connector	Polarographic or optical (option) VP				
pH sensor connector	Gel-filled VP				
Temperature sensor connector	Pt100 with pocket M12				
pH Redox sensor (option)	Gel-filled VP				
Turbidity sensor (option)	Single Channel NIR Absorption Probe				
Pressure transmitter (option)	Piezoresistive M12				
Material (product-wetted parts)	Borosilicate glass Stainless steel AISI 316L/EPDM				

Utility consumption

Power consumption max. DCU Tower Supply Tower	200 2000 [W]
Water consumption max. per Supply Tower	5 L/min
Gas consumption max. per Supply Tower 0.5 L 1 L 2 L 5 L 10 L	1 1.6 4.2 13 20 [L/min]

Utility requirements | Housing connection

Power supply DCU Tower Supply Tower	110 – 230V/6A GFIC: 32mA 120V/15A or 230 V/10A GFIC: 32 mA
Gasses	Controlled @ 1.5 barg (22 psig); dry, particle and oil-free hose connector OD 6 mm
Water	Controlled @ 2–4 barg (29–58 psig) hose connector OD 10 mm
Drain	Gravity drain with zero backpressure required hose connector OD 10 mm
Regulatory compliance	CE (build according to UL&CSA requirements)

Ordering information

Description	BIOSTAT® B-DCU II O ₂ Enrichment				
	0.5 L	1 L	2 L	5 L	10 L
Culture vessel working volume					
Cat. No. 120 VAC					
BIOSTAT® B-DCU II Single	RBD1M5DLOTRDG1	RBD1M01LOTRDG1	RBD1M02LOTRDG1	RBD1M05LOTRDG1	RBD1M1ALOTRDG1
BIOSTAT® B-DCU II Upgrade Kit (Vessel + Supply Tower)	RBD1M5DLOTRDG1E	RBD1M01LOTRDG1E	RBD1M02LOTRDG1E	RBD1M05LOTRDG1E	RBD1M1ALOTRDG1E
Cat. No. 230 VAC					
BIOSTAT® B-DCU II Single	RBD1M5DLOTRDG2	RBD1M01LOTRDG2	RBD1M02LOTRDG2	RBD1M05LOTRDG2	RBD1M1ALOTRDG2
BIOSTAT® B-DCU II Upgrade Kit (Vessel + Supply Tower)	RBD1M5DLOTRDG2E	RBD1M01LOTRDG2E	RBD1M02LOTRDG2E	RBD1M05LOTRDG2E	RBD1M1ALOTRDG2E
Control Unit					
Control Tower					
15" color display with touch screen operation	•				
Automatic pH and pO ₂ calibration routine					
Single probe and group calibration	•				
Control Capabilities Per Vessel					
Temperature, pH, DO (5-stage cascade), Stirrer speed; Foam, Level, Substrate	•				
Vessel Pressure measurement control	○ ○				
Turbidity measurement	○				
Redox measurement	○				
Gravimetric Feed Control	○				
Gravimetric Harvest Control	○				
Extended Password Module	○				
Logbook Module	○				
Supply Tower					
Gas mixing (integrated)	O ₂ -Enrichment				
Rotameter for Sparger [l/min]	•				
Solenoid Valve for O ₂ -Enrichment	•				
Mass Flow Controller					
Total Sparger Flow Air and O ₂	○ ○				
Stirrer Motor	•				
Peristaltic Pumps (integrated)	3				
Feed Pumps speed controlled (integrated external)	○ (up to 3) ○ (up to 2)				
Thermostat system (integrated)	•				
Supervisory Process Control Software					
MFCS/DA for data storage	•				
Culture Vessel					
Jacketed UniVessel®					
Stirrer shaft with Single Mechanical Seal	•				
6-blade disk impeller	2	2	2	2	3
Baffles	-	•	•	•	•
Storage bottles	3	3	3	3	3
Air Inlet and Exhaust filter	2				
Aeration tube with Ring Sparger	•				
Inoculation Addition port	•				
Exhaust Cooler	•				
Addition fitting	2 × 3-way	4-way	4-way	4-way	4-way
Sample- Harvest pipe	•				
Manual sampler	•				
Tray for storage bottles	-	•	•	•	•
pH Electrode, cable	•				
DO Electrode, cable, Clark principle Optical	○ •				
Foam sensor, cable	•				
Level sensor, cable	○				
Temperature sensor Pt 100	•				
Turbidity sensor	○				
pH Redox sensor	○				
Pressure sensor control	○ ○				
Exhaust CO ₂	○				
Tubing, O-Ring (spare set)	•				
Foam Disc (mechanical foam destroyer)	-	○	○	○	○

Broad range of accessories available; Configurable and customizable solutions are available outside of this package. Please contact us for further information.

• = standard, ○ = option, - = not available

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius-stedim.com

Sartorius Stedim Systems GmbH
Schwarzenberger Weg 73-79
34212 Melsungen

Phone +49.5661.71.3400
Fax +49.5661.71.3702

www.sartorius-stedim.com

France

Sartorius Stedim Biotech S.A.
ZI Les Paluds
Avenue de Jouques – BP 1051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI Les Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Franzosengraben 12
A-1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Denmark

Sartorius Stedim Nordic A/S
Hoerskaetten 6D, 1.
DK-2630 Taastrup

Phone +45.7023.4400
Fax +45.4630.4030

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.
Edisonbaan 24
3439 MN Nieuwegein

Phone +31.30.6025080
Fax +31.30.6025099

Spain

Sartorius Stedim Spain SA
C/Isabel Colbrand 10-12,
Planta 4, Oficina 121
Polígono Industrial de Fuencarral
28050 Madrid

Phone +34.91.3586102
Fax +34.91.3588804

Switzerland

Sartorius Stedim Switzerland GmbH
Lerzenstrasse 21
8953 Dietikon

Phone +41.44.741.05.00
Fax +41.44.741.05.09

U.K.

Sartorius Stedim UK Limited
Longmead Business Park
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

America

USA

Sartorius Stedim North America Inc.
5 Orville Drive
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Sartorius Stedim SUS Inc.
1910 Mark Court
Concord, CA 94520

Phone +1.925.689.6650
Toll Free +1.800.914.6644
Fax +1.925.689.6988

Sartorius Stedim Systems Inc.

201 South Ingram Mill Road
Springfield, MO 65802

Phone +1.417.873.9636
Fax +1.417.873.9275

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
Av. Dom Pedro I, 241
Bairro Vila Pires
Santo André
São Paulo
Cep 09110-001

Phone +55.11.4451.6226
Fax +55.11.4451.4369

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente No. 149
Ciudad Satélite
53100 Naucalpan, Estado de México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

China

Sartorius Stedim Beijing
Representative Office
No. 33, Yu'an Road,
Airport Industrial Zone B, Shunyi District
Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim Shanghai

Representative Office
Room 618, Tower 1, German Centre,
Shanghai, PRC., 201203

Phone +86.21.28986393
Fax +86.21.28986392.11

Sartorius Stedim Guangzhou Office

Room 704, Broadway Plaza,
No. 233-234 Dong Feng West Road
Guangzhou 510180

Phone +86.20.8351.7921
Fax +86.20.8351.7931

India

Sartorius Stedim India Pvt. Ltd.
10, 6th Main, 3rd Phase Peenya
KIADB Industrial Area
Bangalore – 560 058

Phone +91.80.2839.1963|0461
Fax +91.80.2839.8262

Japan

Sartorius Stedim Japan K.K.
KY Building, 8-11
Kita Shinagawa 1-chome
Shinagawa-ku
Tokyo 140-0001

Phone +81.3.3740.5407
Fax +81.3.3740.5406

Malaysia

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

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10, Science Park Road, The Alpha
#02-25, Singapore Science Park 2
Singapore 117684

Phone +65.6872.3966
Fax +65.6778.2494

Australia

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

Phone +61.3.8762.1800
Fax +61.3.8762.1828