



BIOSTAT® B-DCU II

Advanced Additive Flow



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 modelling of various culture processes, the BIOSTAT® B-DCU II provides a new level of power and flexibility. The BIOSTAT® B-DCU II Advanced Additive Flow packages are specially configured for cell culture. The integrated automatically-controlled gas mixing system for up to 6 gas flow path provides Sparger and Overlay or optional individual gas out lets. Each gas has its own rotameter for individual flow rate adjustment.

Control Tower

- Graphical user interface with color display and touch screen operation
- Measurement and control for Temperature, pH, DO, agitation and Foam
- User configurable 5-stage DO cascade
- Up to 4 feed controller per vessel
- Automatic gas composition controlled by pH & pO₂ controller
- Level control via Level probe balance
- Totalizers with digital calibration for valves and pumps
- In-process pH-recalibration
- Trend display for up to 8 process values
- Up to 4 balance connections per vessel
- Culture vessel pressure measurement and control

Supply Tower

Advanced Additive Flow Gassing System

- Sparger and Overlay or optional individual gas out lets
- Gasmixing of Air, O₂, N₂, CO₂ for Sparger gassing via solenoid valves
- Air for Overlay gassing
- Additional optional gas flow path
- Optional gas switch Sparger to Overlay
- Up to 6 optional mass flow controller

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
- 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
- 3-blade segment impeller
- Aeration tube with micro Sparger, Overlay aeration fitting, sterile filters and exhaust cooler
- Manual sampler with sampling pipe
- Removable addition bottle support
- 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 BioPAT® DCU control system with 15" TFT color touch screen
- 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 6 Rotameter and Mass Flow Controller
- Culture vessel pressure control
- Easy on-site Supply Tower | Culture vessel upgrade
- Improved connectivity of utilities and probes
- Easy upgrade of cell culture packages for multi-purpose use
- 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 | tbid 20–2000 20–2000 20–1500 20–800 [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 "Advanced Additive Flow" with up to 6 gas flow paths

| | |
|---------------|-------------------|
| Outlet design | Hose tube OD 6 mm |
|---------------|-------------------|

Flowmeter

| | | | | | |
|---|--|-----------|----------|----------|-----------------|
| | Air calibrated @ 1.21 bara 20°C | | | | |
| | 0.5 L | 1 L | 2 L | 5 L | 10 L |
| Gas flow range "Sparger" for Air & N ₂ | 0.53–53 | 0.53–53 | 13–133 | 13–133 | 26–266 [mL/min] |
| Gas flow range "Sparger" for O ₂ & CO ₂ | 1.5–15 | 1.5–15 | 3.3–33 | 3.3–33 | 13–133 [mL/min] |
| Gas flow range "Overlay" for Air | 0.11–1.05 | 0.11–1.05 | 0.16–1.6 | 0.42–4.2 | 1.3–13 [L/min] |
| Additional gas flow path to "Sparger" or "Overlay" | Various flow ranges available from 1.5–15 mL/min to 2.3–23 L/min | | | | |
| Accuracy | +/- 2% FS | | | | |

Thermal Mass Flow Controller (option)

| | | | | | |
|---|--|----------|----------|--------|----------------|
| | Calibrated to specific gas | | | | |
| | 0.5 L | 1 L | 2 L | 5 L | 10 L |
| Gas flow range "Sparger" for Air & N ₂ | 0.5–50 | 0.5–50 | 3–150 | 3–150 | 6–300 [mL/min] |
| Gas flow range "Sparger" for O ₂ & CO ₂ | 0.15–15 | 0.15–15 | 0.6–30 | 0.6–30 | 3–150 [mL/min] |
| Gas flow range "Overlay" for Air | 0.02–1.0 | 0.02–1.0 | 0.03–1.5 | 0.1–5 | 0.2–10 [L/min] |
| Additional gas flow path to "Sparger" or "Overlay" | Various flow ranges available from 1.5–15 mL/min to 0.4–20 L/min | | | | |
| Accuracy | +/- 1% FS | | | | |

Agitation

| | |
|--|-----------------------------|
| Motor performance torque 0.5 L 1 L, 2 L, 5 L, 10 L | 75/0.28 200/1.03 [W]/[Nm] |
|--|-----------------------------|

Integrated pumps

| | |
|--|---|
| Pump head | on off, pulse-width modulated controlled |
| Rotation speed 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 |
| | 0.4–4.4 1.6–16.2 [ml/min] |

Integrated feed pump (option)

| | |
|--|---|
| Pump head | Watson Marlow 102R, for tubing with 1.6 mm wall thickness |
| Rotation speed speed | 1–10 5–50 [rpm] |
| Flow rate tube dependent (bore×wall tubing) 1.6×1.6/3.2×1.6 [mm] | 0.2–2/0.8–8 1.1–11/4–40 [ml/min] |

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 |
| | Quick couplings Quick couplings |

External connections

| | |
|--|-----------------|
| Balance standard option/interface/connector | 2 4/RS232/M12 |
| CO ₂ exhaust analyzer interface/connector | 1/RS232/M12 |
| Feed pump connection qty./interface/connector | 2/0–10 V/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 | 0.15–0.5 L | 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 Power Supply | 5 L/min |
| Gas consumption max. per Power Supply 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 230V/10A GFIC: 32mA |
| 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 Advanced Additive Flow | | | | |
|--|---|-----------------|-----------------|-----------------|-----------------|
| | 0.5 L | 1 L | 2 L | 5 L | 10 L |
| Culture vessel working volume | | | | | |
| Cat. No. 120 VAC | | | | | |
| BIOSTAT® B-DCU II Single | RBD1C5DLATSDG1 | RBD1C01LATSDG1 | RBD1C02LATSDG1 | RBD1C05LATSDG1 | RBD1C1ALATSDG1 |
| BIOSTAT® B-DCU II Upgrade Kit (Vessel + Supply Tower) | RBD1C5DLATSDG1E | RBD1C01LATSDG1E | RBD1C02LATSDG1E | RBD1C05LATSDG1E | RBD1C1ALATSDG1E |
| Cat. No. 230 VAC | | | | | |
| BIOSTAT® B-DCU II Single | RBD1C5DLATSDG2 | RBD1C01LATSDG2 | RBD1C02LATSDG2 | RBD1C05LATSDG2 | RBD1C1ALATSDG2 |
| BIOSTAT® B-DCU II Upgrade Kit (Vessel + Supply Tower) | RBD1C5DLATSDG2E | RBD1C01LATSDG2E | RBD1C02LATSDG2E | RBD1C05LATSDG2E | RBD1C1ALOTSDG2E |
| 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, Substrate | • | | | | |
| Vessel Pressure measurement control | ○ ○ | | | | |
| Turbidity measurement | ○ | | | | |
| Gravimetric Feed Control | ○ | | | | |
| Gravimetric Harvest Control | ○ | | | | |
| Extended Password Module | ○ | | | | |
| Logbook Module | ○ | | | | |
| Supply Tower | | | | | |
| Gas mixing (integrated) | Advanced Additive Flow | | | | |
| Gas outlets | | | | | |
| Sparger & Overlay Individual | • ○ | | | | |
| Rotameter Sparger | • Air, O ₂ , N ₂ , CO ₂ | | | | |
| Rotameter Overlay Additional gas | • Air ○ O ₂ or CO ₂ or N ₂ | | | | |
| Solenoid valves for gas mixing of Air, O ₂ , N ₂ , CO ₂ | • | | | | |
| Gas switch Sparger to Overlay | ○ | | | | |
| Mass Flow Controller | ○ (up to 6) | | | | |
| Stirrer Motor | • | | | | |
| Peristaltic Pumps (integrated) | 3 | | | | |
| Feed Pumps speed controlled (integrated external) | ○ (up to 3) ○ (up to 2) | | | | |
| Thermostat system (integrated) | | | | | |
| Jacketed vessel Single wall vessel | • ○ | | | | |
| Supervisory Process Control Software | | | | | |
| MFCS/DA for data storage | • | | | | |
| Culture Vessel | | | | | |
| Jacketed vessel Single wall vessel | • - | • ○ | • ○ | • ○ | • ○ |
| Stirrer shaft with Single Mechanical Seal | | | | | |
| Direct coupling Magnetic coupling | • - | • ○ | • ○ | • ○ | • ○ |
| Magnetic drive | - | - | ○ | ○ | ○ |
| 3-blade segment impeller | 1 | | | | |
| Storage bottles | 3 | 3 | 3 | 3 | 3 |
| Air Inlet and Exhaust filter | 3 | | | | |
| Aeration tube with μ-Sparger Overlay port | • | | | | |
| 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 | ○ | | | | |
| Exhaust CO ₂ sensor | ○ | | | | |
| Pressure sensor control | ○ ○ | | | | |
| Tubing, O-Ring (spare set) | • | | | | |

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