

# Software Features

## BioPAT® MFCS

### Benefits

- Scalable software for almost all bioprocess applications
- Central platform for real-time data and historians of process, analytical and sampling data
- Automated batch execution with the ANSI/ISA-88 compliant recipe control module
- Supports all requirements to archive GMP and FDA 21 CFR Part 11/EU Annex 11 compliance
- Fully user configurable and upgradeable with specific modules
- Installation, configuration, validation and engineering services

### Product Information

BioPAT® MFCS software enables you to incorporate a new standard in bioprocess data management and automation.

Its reliable data acquisition, efficient trend monitoring, and advanced recipe control make it an ideal tool for all upstream and downstream processes, no matter if you prefer single-use or reusable systems. The new BioPAT® MFCS is your solution for robust and reproducible processes in R&D and GMP environments – all backed by our expertise since 1986.



# BioPAT® MFCS/win Support and Lifecycle

Main Support End: 13<sup>th</sup> October 2024

Hotline support, delivery of patches for technical or security issues

Extended Support End: 14<sup>th</sup> October 2026

End of Hotline Support

## BioPAT® MFCS Software Features

Feature	BioPAT® MFCS	BioPAT® MFCS/win 3.1
Data acquisition and storage	■	■
Supervisory control	■	■
Process value alarms	■	■
Process value simulation	■	■
Configuration management	■	■
Sample data management	□	■
Calculations	□	■
Advanced charting	■	■
Process Unit Display	■	■
<b>Recipes</b>		
Recipe control according to ISA-88.01	□	□
Perfusion Recipe	□	□
Customized Recipes	□	□
<b>Networking</b>		
Multi User Client	□ (limited to 30 clients)	□ (limited to 10 clients)
<b>GMP &amp; User Management</b>		
21 CFR Part 11 compliance	□	□
User management	□	□
<b>Others</b>		
Replication manager	■	□
Integrated Multivariate Data Analysis (MVDA)	■ (SIM API)	□
Virtualized Environment	■	■
Feature	BioPAT® MFCS	BioPAT® MFCS/win 3.1
<b>Connectivity</b>		
OPC DA Client	□ (R/W; Classic & UA)	□ (R/W; Classic)
OPC Classic DA server	■ (requires UA gateway)	□
OPC UA server (DA&AC)	□	-
OPC Classic AE server	■ (requires UA gateway)	□
Remote Alarming Package	□	□

<b>General</b>		
Number of supported process units	up to 32	up to 16
Minimum data export interval	0.5 s	1 s
Supported database systems	Microsoft® SQL 2019 and 2022	Microsoft® Access
User interface languages	EN, DE, ZH, FR, ES, IT	EN
<b>Training &amp; Support</b>		
Training	<input type="checkbox"/>	<input type="checkbox"/>
Validation support	<input type="checkbox"/>	<input type="checkbox"/>
Service and Support Agreement	■ (free of charge in the first year)	<input type="checkbox"/>
<b>Supported Operating Systems</b>		
Windows® 10 <sup>1</sup>	32 bit   64 bit	64 bit
Windows® 11 <sup>1</sup>	32 bit   64 bit	-
Windows® Server 2016	64 bit	64 bit (only up to version Windows® 10 Pro 64 bit)
Windows® Server 2019	64 bit	64 bit
Windows® Server 2022	64 bit	-
■ included <input type="checkbox"/> optional		

Microsoft® and Windows® are registered trademarks of Microsoft Corporation, USA

<sup>1</sup> No support for LTSC/LTSB, Home version and versions no longer supported by Microsoft®

## Minimum System Requirements for BioPAT® MFCS

### Server & Client Stand-alone Installation

- 8-core processor 3 GHz
- 32 GB memory (RAM)
- 500 GB hard disk space
- DirectX, 1GB RAM graphic card
- LAN adapter, RJ-45, 1 Gbit
- 15.6" monitor (1366 × 768)

### Supported Virtual Environments

- VMware® Workstation
- Hyper-V®

### Multi-User Installation

#### Client

- 6-core 3 GHz processor
- 16 GB memory (RAM)
- 50 GB hard disk space
- DirectX, 1GB RAM graphic card
- LAN adapter, RJ-45, 1 Gbit
- 15.6" monitor (1366 × 768)

#### Server


- 6-core 3 GHz processor
- 32 GB memory (RAM)
- 500 GB hard disk space
- LAN adapter, RJ-45, 1 Gbit
- DirectX, 1GB RAM graphic card

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

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