

BioPAT® MFCS | Win Manager Training

Become an Expert and use
the full potential of your
BioPAT® MFCS Software



Training Content

Overview about BioPAT® MFCS | win and its basic functions:

- Operator Service Program, Data Management, Configuration options
- Configuration of Process Units and Control Units
- PID Control and Applications
- User Management, 21 CFR Part 11 and Audit Trail
- Configuration Reports
- Update Configuration and Backup Copy
- Maintenance, System Recovery and System Utilities
- Processes automation with S88 Recipes
- Installation and Uninstallation of BioPAT® MFCS | win Server | Workstation

- BioPAT® MFCS | win Network Structure and Settings
- Security Settings- Help-Documentation

Training Setup


- Group size: Up to 8 participants
- Where: At customer side
- Course duration: 3 days

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