

BioPAT® MFCS | Win Basic User Training

Discover the possibilities with
the BioPAT® MFCS Software



Training Content

- Introduction into SCADA-System BioPAT® MFCS | win
- Shell and Service Control Program
- Data storing and Batch Management
- Process Visualization with Operator Service
- Batch Start, Batch Setup and Initialize defined Batch
- Process automation with S88 Recipes (Using and Modification)
- Off-line Data Management
- Evaluation by Plotting
- Reporting of Batches and Data export
- 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