

BioPAT® MFCS | Win Trouble Shooting Training

Help yourself and become an Expert on the functionalities of your BioPAT® MFCS Software



Training Content

- Crossflow filtration, filtration parameters
- Transmembrane pressure, Reynolds-number, yield
- Introduction into SCADA-System BioPAT® MFCS | win
- Shell, Service Control Program
- Data storing and batch management
- Process visualization with Operator Service Program
- Off-line data management
- Evaluation by plotting- Reporting of batches and data export
- Configuration of variables, controllers and calculations
- DCU sequences for filtration processes
- Configuration of equipment phases
- Update configuration und backup copy
- Automatization of filtration processes with S88 recipes

Training Setup


- Group size: Up to 8 participants
- Where: At customer side
- Course duration: 2 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