

BioPAT® MFCS|4 Basic User Training

Discover the possibilities
within the BioPAT® MFCS
Software



Training Content

- Getting started with BioPAT® MFCS 4
- Configuration of devices and units
- Process data acquisition and visualization
- Creation of new variables, online and offline calculations
- Visualization of calculated values and offline variables
- Charting of live and historical data
- Introduction into process automation with the Recipe Control Module
- Creation and modification of ISA-88 compliant recipes
- Operations and Phases
- Branches and conditions
- Event and time-based transitions
- Execution of recipe controlled batch processes

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