

# BioPAT® MFCS | Win Advanced User Training

Get the best out of your  
BioPAT® MFCS Software



## Training Content

- Advanced Training BioPAT® MFCS | win
- Integration, Differentiation
- MIN, MAX Determination
- Moving Average, Timers
- If then Else Structures
- OUR, CER, RQ Calculation
- Exponential Feed
- PID Controllers and MFCS PID Control
- pO2 Cascade Control (DO Control)
- Gravimetric Flow Control Fedbatch Fermentation with Gravimetric Flow Controllers

- Continuous Fermentation with Gravimetric Flow Controllers
- DO-Stat | pH-Stat
- kLa | Determination of kLa

## 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](http://www.sartorius.com)