

# BioPAT® MFCS | 4 Trouble Shooting Training

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



## Training Content

### **The knowledge transfer training that enables the Operator or Administrator to:**

- Use the MFCS Shell and Service Control Program
- Create an MFCS Database Structure
- Use Convert Tools (Import Batches, Update Databases)
- Work with the Diagnostic Tools (System Diagnostic, Maintenance Tool)
- Make use of the Repair Tools (Re-installation Dongle driver, Repair of Databases)
- Deploy the Access 2000 Runtime installation (ART\_2000)
- Change Windows Settings and Network Installation
- Carry out Database Backup and System Recovery
- Maintain the PC hardware

- Maintain the operating system (Hard disk defragmentation)
- Make use of the Archiving function of the batch database
- Move batches to archive location
- Enable Maintenance function for "Replicated Batches", if Replication Manager is installed

## Training Setup


- Group size: Up to 8 participants
- Where: At customer side
- Course duration: 1 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)