## SARTURIUS

# BioPAT® MFCS | 4 Remote Training

Discover the possibilities within the BioPAT® MFCS Software.



## **Training Content**

#### MFCS Overview

- Database Management
- Server-Client
- Network
- Help Documents

#### Administration

- Configuration
- Calculations
- Recipe S88 Module
- Back Up

#### Monitoring

- Batch Start and Stop
- Trend-, Unit-, Alarm- and Recipe Display

### Analysis:

- Charting of historian Batches
- Sample Data Management
- Data Export and Report

## **Training Setup**

Group size: Up to 8 participants

• Where: Remote e.g. via MS Teams

Duration: 8 hours max

#### Germany

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen Phone +49 551308 0

For further contacts, visit www.sartorius.com