

Service & Support Package

BioPAT® MFCS 4

Benefits

- Efficient technical support and troubleshooting
- Quick response
- Reliable system performance



Product Information

Sartorius prides itself on the quality and reliability of our BioPAT® MFCS software. Like all professional software, dedicated services are required to ensure the most efficient performance over the long term. This is why we are pleased to offer you our comprehensive, worry-free Service & Support Package. Maximize the performance of your BioPAT® MFCS software and contact our Service Team today!

Your one-year Service & Support Package includes all of the following:

- Direct access to our Service Team by e-mail with timely response by e-mail or phone
- Remote desktop assistance
- Downloads of product updates
- Downloads of security updates
- Access to free training webinars
- Newsletters with information updates on upcoming BioPAT® MFCS modules
- Options to extend your Service & Support Package

Your direct link to our Service Team: mfcs.service@sartorius.com

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen, Germany Phone +49.551.308.0 Fax +49.551.308.3289 www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178 Argentina +54.11.4721.0505 Brazil +55.11.4362.8900 Mexico +52.55.5562.1102 UK +44.1372.737159 France +33.442.845600 Italy +39.055.63.40.41 Spain +34.913.586.098 Russian Federation +7.812.327.53.27 Japan +81.3.4331.4300 China +86.21.6878.2300



