

Interface Consulting

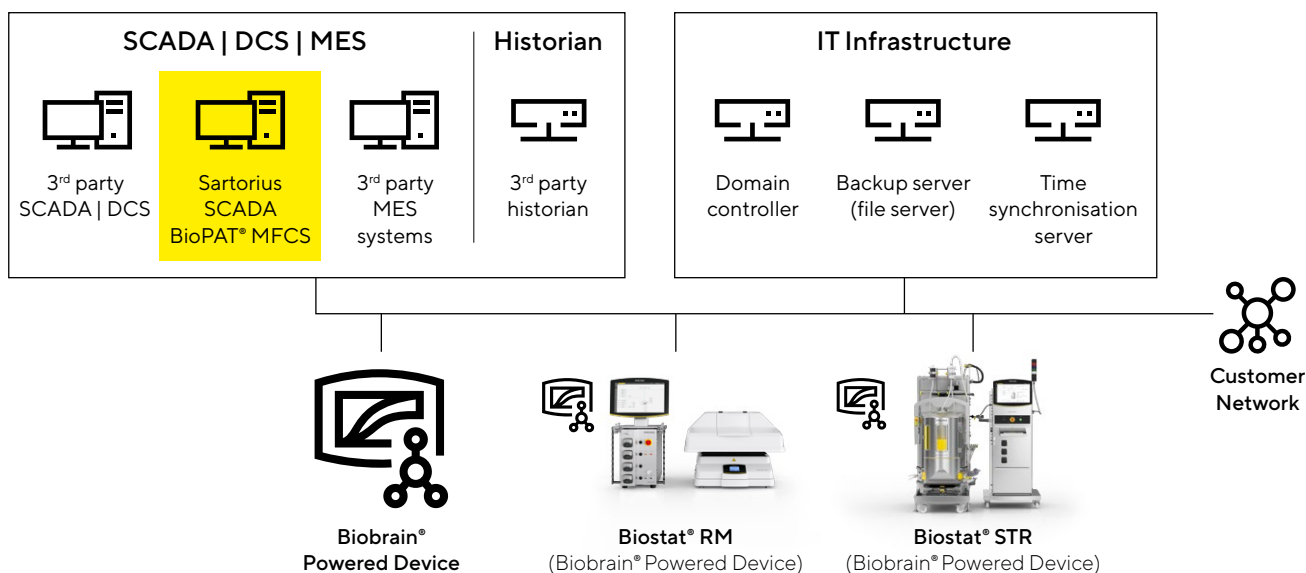
Support and Expertise to Enhance Your Success

Integrating Biobrain®-powered products and | or BioPAT® MFCS software into your higher level systems (for example, IT infrastructure or DCS) can be complex and time-consuming. Our Interface Consulting Services provide expert guidance, specialized knowledge transfer, and intuitive integration documentation to overcome implementation challenges to mitigate risk, reduce effort, increase productivity, and shorten timelines.

- Virtual consulting time with experts for training, specialized knowledge transfer, support, and Q&A
- Access to intuitive integration documentation, including pre-verified, step-by-step configuration instructions for typical integration scenarios
- Demonstration of typical integration scenarios in a virtual environment
- Support for higher level systems, including Siemens PCS7®, Emerson DeltaV®, Rockwell PlantPAx® DCS, and BioPAT® MFCS



Order Now



Consulting Service for Integrating BioPAT® MFCS Into Facility Environment

Attribute	BioPAT® MFCS
Remote consulting time with dedicated engineer	8 hours
Domain integration and network setup	✓
Instrument and software name definition setup	✓
Security (antivirus, communication)	✓
Backup & restore data and system	✓
Number of Instruments that can be connected per BioPAT® MFCS	1 - 32
Instruments supported per consulting engagement (order)*	Up to 32
MFCS systems supported per consulting engagement (order)*	5

Ordering Information

Order Number:
ZB-MB-HA-0044

Sartorius' DCS Partners



Consulting Service for Integrating Biobrain® Powered Products

Attribute	Basic Small	Basic Medium	Basic Large	Advanced Small	Advanced Medium	Advanced Large
	1-3 Instruments	4-12 Instruments	13-24 Instruments	1-3 Instruments	4-12 Instruments	13-24 Instruments
Description	Provides consulting and documentation for network integration			Provides consulting and documentation for network integration and DCS integration via OPC UA		
Remote consulting time with dedicated engineer	8 hours	16 hours	24 hours	16 hours	24 hours	36 hours
Domain integration and network setup	✓	✓	✓	✓	✓	✓
Instrument and software name definition setup	✓	✓	✓	✓	✓	✓
Security (antivirus, communication)	✓	✓	✓	✓	✓	✓
Backup & restore data and system	✓	✓	✓	✓	✓	✓
Data Acquisition from to Biobrain® into DCS for visualization, supervisory control, Historization, alarms and events, and alarms acknowledgement				✓	✓	✓
Recipe execution in DCS via phase-level integration of Biobrain® powered product equipment phases				✓	✓	✓
Recipe level integration in DCS by triggering recipes in the Biobrain® powered product via DCS****				✓	✓	✓
Siemens PCS7 DCS**	✓	✓	✓	✓	✓	✓
Emerson DeltaV DCS****	✓	✓	✓	✓	✓	✓
Rockwell PlantPAX® DCS**	Please Contact your Sales Representative					
Biobrain®-powered products (e.g., Biostat® RM, Biostat STR® GEN 3)	✓	✓	✓	✓	✓	✓

*Mix and match up to 32 instruments distributed in a maximum of 5 MFCS systems

**Consulting service for one DCS platform per engagement

***Biobrain® version 8 and higher

****Biobrain® version 8 and higher. Not available for DeltaV

*****Currently, demonstration of typical integration scenarios in a virtual environment available for Siemens PCS7 only