



BioPAT® MFCS/win

Recipe Control (S88) Module

Product Information

The time has come for robust processes – the BioPAT® MFCS/win Recipe Control (S88) module:

- Improved batch-to-batch consistency
- Enhanced process understanding
- Standardized automation
- Decreased risk of errors
- Simplified tech-transfer
- Minimized number of rejected lots
- Free up operator time

Introduction

BioPAT® MFCS/win is the world standard software for supervisory bioprocess control and data acquisition. Based on over 25 years of experience and more than 3,000 installations, the latest release of BioPAT® MFCS/win provides various solutions to meet your specific requirements.

Besides the core functionality of a true SCADA system for reusable and single-use bioprocess applications, BioPAT® MFCS/win offers a number of additional modules for adaptation to individual requirements. The unique Recipe Control (S88) Module allows bioprocess operators to realize automation strategies for batch, fed-batch and continuous processes in-line with Quality by Design (QbD) principles.

The American ANSI/ISA-88.01 standard and its international equivalent IEC-65112 are industrial standards for the control of all kinds of batch processes. The BioPAT® MFCS/win Recipe Control (S88) Module follows these standards and contains all functions and models that are of use for biopharmaceutical up- and downstream processes.

Minimize run-to-run variability

The BioPAT® MFCS/win software package as well as the Recipe Control (S88) Module support industrial users as well as researchers in universities and institutes. Whereas the industrial user at the



pilot or production plant may need a well-structured, validatable system, the researcher needs a very flexible and open system. The Recipe Control (S88) Module considers both aspects.

In order to achieve a reproducible and fully automated batch process, BioPAT® MFCS/win recipes are designed to trigger different actions automatically by events. Initialization of substrate feed control after automatic detection of the end of a batch or to automatic start of harvesting after a specific cell concentration has been reached are only two examples.

The S88 recipe structure adapts perfectly to constantly evolving recipes in the pilot plant development area. As it provides a common language for process design and specification, an early recipe development enables the implementation of an operator independent and robust process control strategy across development and commercial manufacturing.

The advanced use of Process Analytical Technologies (PAT), whether in development to gain process understanding or manufacturing to monitor Critical Process Parameters (CPPs), enables the implementation of enhanced control strategies. With its conditional logic and the direct translation of biopharmaceutical unit operations into S88 phases the Recipe Control Module supports your automation strategy to reduce variability and thus increase batch-to-batch consistency.



Select the pointer to move phases with drag & drop



Switch on variables and data storage, set alarm limits, change controller properties, start calculations, etc.



Run a time profile and change variable value, controller setpoint or output



Define a controller setpoint and verify that value stays within the defined deadband for a certain time



Start any program (*.exe) and optionally hand over parameters



Start a DCU Sequence which is defined in the configuration of the digital control unit



Enter a prompt message, which will pop-up during recipe execution



Link 2 process units which depend on each other



Features

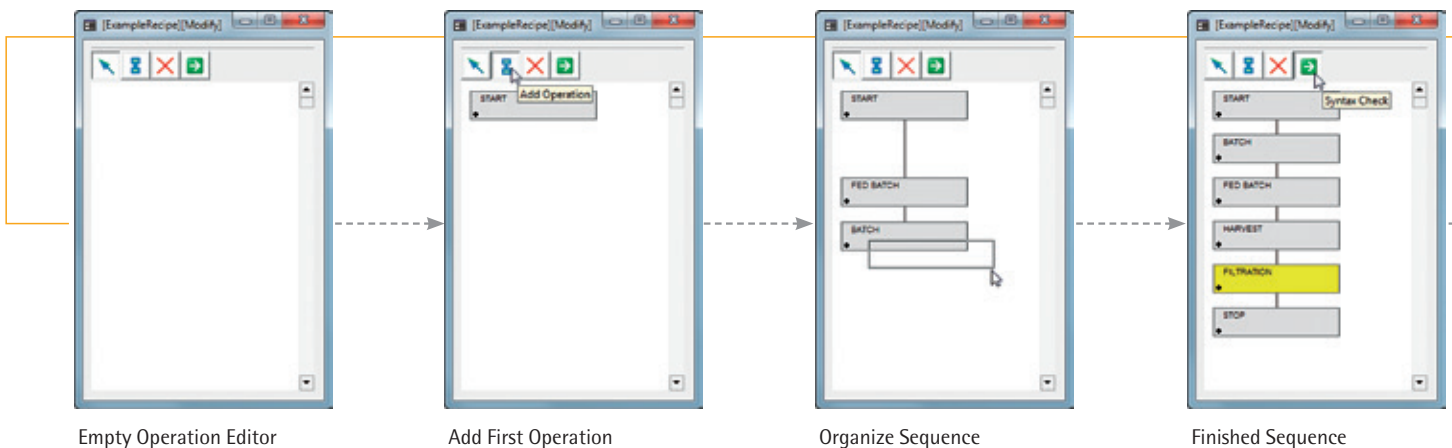
- Graphical Configuration and Execution
- Fully mouse-driven recipe editor
- Drag & drop for operations and phases
- Comprehensive toolbar with pre-defined phases
- Color coded view of current process state
- Operation and Phase library
- Automatic Syntax Check
- Automatic Recipe Version Control
- Possible restriction to equipment-classes
- Review & approve functionality

Easy Graphical Programming

The new recipe editor permits graphics-based process automation with loops and jumps, facilitating application-oriented automation in compliance with industry standards. Semi or fully-automatic operations and pre-defined phases with state or time-dependent transitions enable an organized and structured batch processing as well as flexible manufacturing.

Efficient Workflow Optimization with DoE

A key benefit of applying QbD is the ability to identify critical quality attributes and process parameters. The correlation can be made in a cost effective and statistically significant manner by using Design of Experiments (DoE). Based on your experimental objective the unique and additionally available BioPAT® MFCS/win DoE Module generates an appropriate design and automatically configures a S88 recipe.

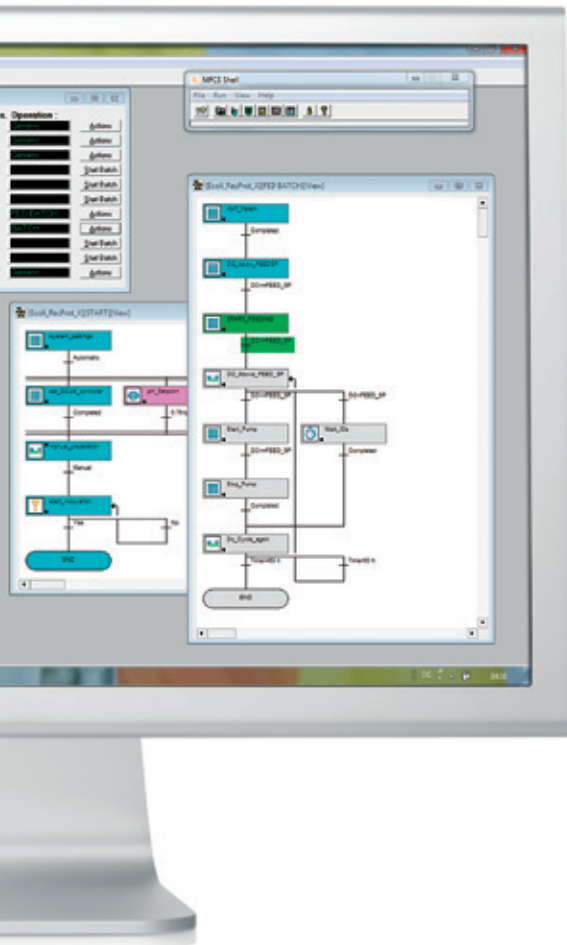










Empty Operation Editor

Add First Operation

Organize Sequence

Finished Sequence



-  Empty phase without any function, e.g. as input phase for jumps
-  Timer, which will be active until a certain day time or for a given duration
-  Every operation must have exactly one End Phase at the last position
-  Create transitions and connect up to 4 phases, where only one single branch becomes active
-  Create multipliers that connect up to 10 phases, where all parallel branches start simultaneously
-  Use jumps to repeat an entire or only certain parts of an operation
-  Delete transitions, multipliers, branches, phases, etc.
-  Mandatory plausability check for errors in order to use a master recipe as control recipe



Faster Tech-Transfer to Manufacturing

The Recipe Control (S88) Module gives you the possibility to center your process development activities around a recipe structure with a common terminology. As recipes scale-up and become more detailed, S88 is ideal for handling the various stages from research to development to commercial manufacturing. Standardized approaches and straightforward mechanism across the organization will finally lead to a simplified technology transfer.

Work in-line with QbD Principles

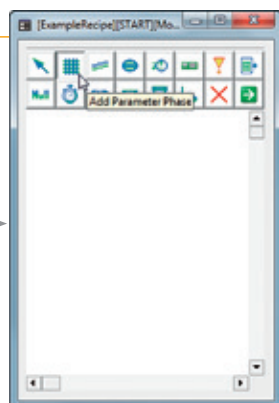
One of the greatest benefits of implementing QbD is the increased regulatory flexibility as well as a faster regulatory approval of new product applications and process changes. BioPAT® MFCS/win Recipes support the realization of advanced scientific concepts and the implementation of enhanced control strategies with global regulatory acceptance.

Professional Service & Support

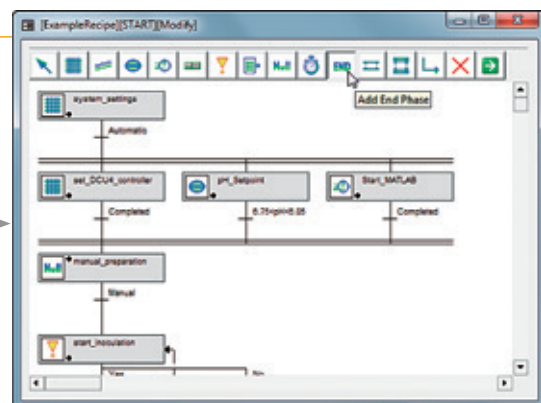
Our team of experienced service engineers has been involved in a number of successful projects world-wide. Expertise gained through long term cooperation with pharmaceutical and biotech companies has been used to develop and refine a comprehensive range of services including computer system validation (CSV) as well as S88 recipe programming based on your specific requirements. In addition the BioPAT® MFCS/win Manager or Application Training provide an excellent basis to get started creating your own recipes.



Modify Operation "Start"



Empty Phase Editor



Add Phases, Timers, Promts, Jumps, Conditions

Applications

BioPAT® MFCS/win provides connectivity to both upstream and downstream devices, so that the Recipe Control (S88) Module is widely used to establish a sophisticated control strategy across both areas and even link systems for integrated bioprocessing to optimize yield and cost-effectiveness of production.

The first common application relates to event detection, which means the identification of process states based on online measurements.

- Automatic batch end detection, e.g. based on DO/offgas measurements
- Initialization of feed | induction, e.g. based on biomass measurement
- Just-in-time harvest, e.g. based on (by-) product measurement
- Automatic stop of diafiltration, e.g. based on conductivity measurement

Furthermore the implementation of advanced feeding strategies for fed-batch bioreactors can be realized.

- Linear, exponential or pulse-wise feeding of growth and production culture media
- DO-stat or adaptive control techniques
- Open-loop or closed-loop control of cell specific growth rate μ

Finally recurrent events can be implemented based on loops and jumps.

- Intermittent medium addition or harvesting
- Activation or deactivation of controllers based on user-defined intervals
- Optimization of cross flow performance by periodic adjustment of differential pressure



Gold Application Development

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

Specifications subject to change without notice.
Copyright Sartorius Stedim Biotech GmbH. Printed
in the EU on paper bleached without chlorine.
Publication No.: SLL2008-e170403
Order No.: 85032-544-99
Ver. 04 | 2017