



Rapid and scalable feed preparation with powdered soy hydrolysates or yeastolates and single-use magnetic mixing systems



Application
Note

#18

#19

#20

#21

#22

Executive summary

This application note presents a new approach to increase the speed and consistency of feed preparation steps in bio manufacturing. The method combines ready to use feed formulations from SAFC[®], with scalable, high efficiency, single-use mixing systems provided by Sartorius Stedim Biotech.

Two examples of feed preparation steps, 50X BD Yeastolate and 50X SE50MAF-UF Soy Hydrolysate, are presented for different volumes. The contained transfer of powdered feed formulations into the high torque single-use mixing system enables a rapid dissolution and dispersion of the feed powders in liquid for volumes of 50 L and 200 L. The performances of the single-use mixing system are characterized with quantitative (conductivity measurement) and qualitative (visual inspection) techniques. The proposed method and system provide seamless scale-up and consistent rapid feed mixing for process development and GMP manufacturing.

Introduction

This application study presents the performances of a fully single-use mixing solution for the large scale preparation of two different feeds. The prepared feed are concentrated 50 times and can therefore be used for large volume media preparation using the appropriate dilution. The mixing technology selected for this application is Flexel[®] for Magnetic Mixer¹ with volumes of 50 L and 200L. The magnetic coupling of the impeller with the Magnetic Mixer Drive Unit enables a rotation speed up to 300 rpm, providing a powerful mixing of the feed.

The first feed tested in this study was the 50X BD Yeastolate. This feed is formulated specifically for media used with insect cell culture and cell lines that utilize L-Glutamine. It is generally used at a final formulation of 4 g/L.

The second feed tested in this study was 50X SE50MAF-UF Soy Hydrolysate. This feed is an animal-component free feed utilized in cell culture and fermentation applications as a nitrogen source when non-animal components are required. It is generally used at a final concentration of 5 g/L for cell lines which do not need any glutamine addition like CHO-GS.

Purpose of the application study

The purpose of this application study is to assess the performances of the Flexel[®] for Magnetic Mixer technology to dissolve the two feeds. The mixing times are determined by conductivity and visual inspection of the solution in the Flexel[®] Bag for Magnetic Mixer.

Materials and methods

The list of materials and equipments used for this application is:

1. Standard Flexel[®] Bags for Magnetic Mixer (50 L: FMB114867, 200 L: FMB114893).
2. Powder Transfer Bag System (15 L : FMA114008, 30 L: FMA114009).
3. Palletank for LevMixer[®] | Magnetic Mixer (50 L: FXC110820, 200 L: FXC110821).
4. Magnetic Mixer Drive Unit, 230 V, EU power cord (ref. LT-DU-006-EU).
5. Powder Bag holder (ref. FXA114344).
6. SAFC[®] feed:
 - 50X SE50MAF-UF Soy Hydrolysate (Product number: 58903C/2019219)
 - 50X BD Yeastolate (Product number: 58902C/2025050)
7. Conductivity sensor: WTW InoLab Cond 740i.
8. Floor scale: Sartorius IF S4 1500RR-1.

Method used:

1. The Flexel[®] Bag for Magnetic Mixer is placed into the Palletank with conductivity and temperature sensors.
2. The bag is filled with deionised water to 80% of the nominal volume (water temperature: 20 °C).
3. The mixing speed is turned on and set up to the maximum speed of 300 rpm to optimize powders dispersion.
4. Feed powders are added slowly through the top port to ease the powder incorporation into the water
 - 50X BD Yeastolate
 - 50X SE50MAF-UF Soy Hydrolysate

The powders are incorporated in the Flexel[®] Bag for Magnetic Mixer using either:

- SAFC[®] bucket liner
- or Sartorius Stedim Biotech Powder Transfer Bag for a contained transfer to the mixing bag assembly

The actual concentrations of feed during the mixing phase are:

- 50X BD Yeastolate : 214 g/L (density = 1.068 kg/L)
- 50X SE50MAF-UF Soy Hydrolysate : 257 g/L (density = 1.078 kg/L)

¹ This product uses Pall patented Magnetic Mixer technology. All information on patents can be found at Pall.com/patents.

5. Two mixing times are monitored from the addition of feed powders:

5.1 "mixing time 1" is determined from the conductivity signal as follows:

The "mixing time 1" corresponds to the time when 95% of the final value is reached and when all next measurements stay within a 5% tolerance.

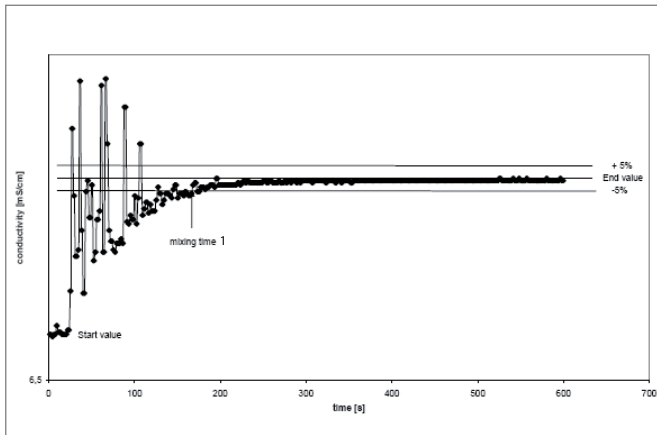


Figure 1: General principle of mixing time determination via conductivity

5.2 "mixing time 2" is determined by a visual inspection. The "mixing time 2" corresponds to the time when all suspended particles are visually dissolved.



200 L Flexel® with Magnetic Mixer Technology



200 L Palletank for Magnetic Mixer equipped with the Powder Transfer Bag System for the mixing trial

Results and discussions

1. Mixing performances results

50X BD Yeastolate and 50X SE50MAF-UF Soy Hydrolysate feed have been prepared in Flexel® Bag for Magnetic Mixer at 50 L and 200 L scales. Mixing time's results are presented for each feed at the different scales.

Mixing time results for 50X BD Yeastolate – 50 L

For 50X BD Yeastolate preparation at 50 L, 10 kg of powder were added in less than 2 minutes in the mixing bag using SAFC® bucket liners. The feed powder was mixed in less than 8 minutes (visual check).

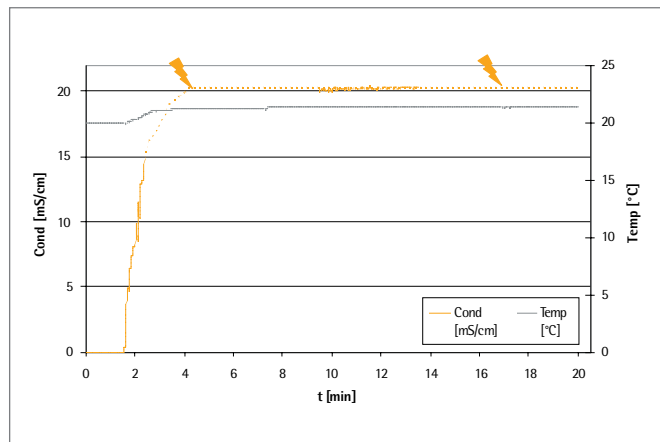


Figure 2: 50X BD Yeastolate preparation in 50 L Flexel® Bag for Magnetic Mixer

Note: conductivity values are not stable at 50 L due to the presence of air bubbles around the conductivity cell (strong vortex). A stable value of the conductivity could be observed only at lower impeller rotation speed.

Mixing time results for 50X BD Yeastolate – 200 L

For the 50X BD Yeastolate preparation at 200 L, 40 kg of powder were added in less than 6 minutes into the mixing bag using two Powder Transfer Bags. The feed powder was mixed in 15 minutes according to conductivity and in 20 minutes according to visual check.

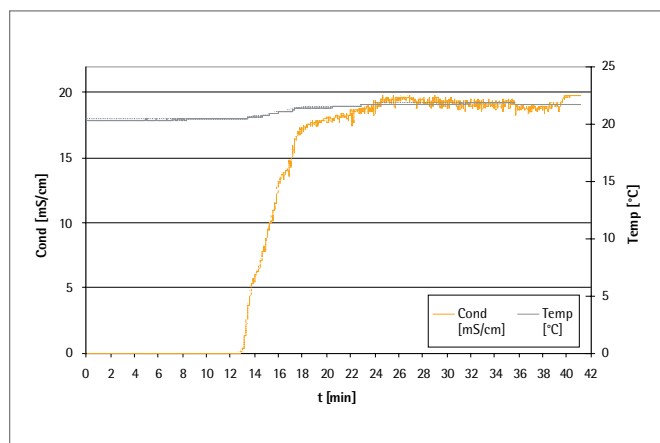


Figure 3: 50X BD Yeastolate preparation in 200 L Flexel® Bag for Magnetic Mixer

Mixing time results for 50X SE50MAF-UF

Soy Hydrolysate – 50 L

For the 50X SE50MAF-UF Soy Hydrolysate preparation at 50 L, 12.5 kg of powder were added in less than 5 minutes in the mixing bag using SAFC® bucket liners. The feed powder was mixed in less than 20 minutes (visual check).

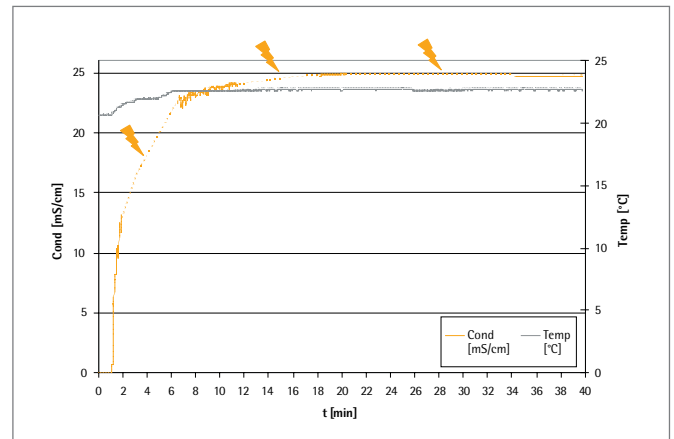


Figure 4: 50X SE50MAF-UF Soy Hydrolysate preparation in 50L Flexel® Bag for Magnetic Mixer

Note: conductivity values are not stable at 50L due to the presence of air bubbles around the conductivity cell (strong vortex). A stable value of the conductivity could be observed only at lower impeller rotation speed.

Mixing time results for 50X SE50MAF-UF

Soy Hydrolysate – 200 L

For the 50X SE50MAF-UF Soy Hydrolysate preparation at 200 L, 50 kg of powder were added in 15 minutes in the mixing bag using five Powder Transfer Bags. The feed powder was mixed in 30 minutes according to conductivity and 47 minutes according to visual check.

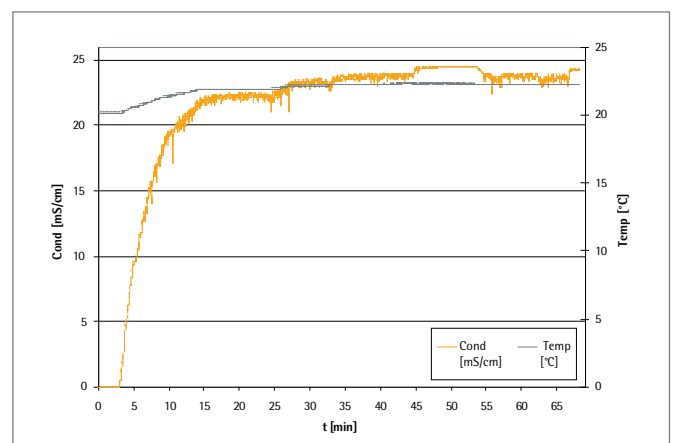


Figure 5: 50X SE50MAF-UF Soy Hydrolysate preparation in 200 L Flexel® Bag for Magnetic Mixer

Mixing performances vs. feed type

The mixing time reported in the table represents only the time to dissolve the feed powders at the different scales. These mixing times include the transfer time of the multiple Sartorius Stedim Biotech Powder Transfer Bags System (for the 200 L scale) or SAFC® bucket liners (for the 50 L).

Volume (L)	Control test	50	200
50X BD Yeastolate	Conductivity Mixing time 1		< 15 min
	Visual inspection Mixing time 2	< 8 min	< 20 min
50X SE50MAF-UF Soy Hydrolysate	Conductivity Mixing time 1		< 30 min
	Visual inspection Mixing time 2	< 20 min	< 47 min

Figure 6: Overview on mixing times for the two powder feed dissolution

General comments

- A rapid dissolution was observed for both 50X SE50MAF-UF Soy Hydrolysate and 50X BD Yeastolate feed due to the strong mixing torque at 300 rpm.
- For the 50 L volume, the vortex at 300 rpm in the 50 L bag volume resulted in the generation of air bubbles that interfered with conductivity measurement with air trapped in the conductivity cell. The conductivity was not unwavering even though mixing was completed. A stable value of the conductivity could be observed only at lower impeller rotation speed. Therefore mixing time cannot be precisely defined using the conductivity signal. Mixing times were determined by visual inspection.
- For the 200 L volume, the conductivity of the solutions reaches a stable value. However, some fine particulates can still be visually observed in the solution. The agitation at 300 rpm was maintained until the particulates became visually totally dissolved. This visual control is facilitated by the large windows of the Palletank.
- A 4x increase in feed preparation volume only leads to a 2x increase in the mixing times.
- The worst case feed preparation condition (200 L 50X SE50MAF-UF Soy Hydrolysate) with mixing of 50 kg of powder in 160 L of water resulted in a mixing time of only 47 min.

Conclusions

- Large volume feed solutions are quick and easy to prepare using the combination of ready to use feed formulations and the high efficiency mixing of the Flexel® with Magnetic Mixer technology.
- Such high concentrated feed preparations are useful in large scale bioproduction. For instance, a 200 L feed preparation at 250 g/L used at 5 g/L in media formulations enables to work with a 10,000 L bioreactor.
- The contained processing conditions with the closed Powder Transfer Bag System docked onto the sterile Flexel® Bag for Magnetic Mixer are favourable to maintain low bioburden and to reduce to the minimum exposure of the operator to chemicals.
- The platform provides a single-use scalable feed preparation capability with a range of Flexel® Bags including volumes of 50 L, 100 L and 200 L.

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany
Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49.551.308.0

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen
Phone +49.5665.407.0

France
Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria
Sartorius Stedim Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium
Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary
Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy
Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands
Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
filtratie.nederland@sartorius-stedim.com

Poland
Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.647.38.40

Russian Federation
LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain
Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland
Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.
Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine
LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA
Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina
Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil
Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico
Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Peru
Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima
Phone +51.1.441 0158

Asia | Pacific

Australia
Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China
Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India
Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan
Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300

Malaysia
Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore
Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea
Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700



▶ www.sartorius-stedim.com