



Rapid and scalable buffer preparation with single-use magnetic mixer



Application
Note

#13

#14

#15

#16

#17

Executive Summary

This application note presents a new approach to increase the speed and consistency of buffer preparation steps in bio manufacturing. The method combines ready to use buffer formulations from SAFC[®], with scalable, high efficiency, single-use mixing systems provided by Sartorius Stedim Biotech. The mixing system is a key component of FlexAct BP, a configurable disposable solution for buffer preparation.

Three examples of large volume buffer preparation steps (Tris | Tris HCl | NaCl, citric acid | sodium citrate | NaCl, PBS), in nominal (x1) or concentrated (x20) forms are presented. The contained transfer of powdered buffer formulations into the high torque single-use mixing system enables a rapid dissolution and dispersion of the buffer powders in liquid for volumes of 50 L to 1,000 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 buffer 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 three different buffers. The mixing technology selected for this application is Flexel[®] for Magnetic Mixer¹ with volumes of 50L, 200L and 1,000L. 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 buffer salts. The three buffer solutions prepared in this application study (Citric acid | sodium citrate, PBS and Tris buffer) are commonly used in biomanufacturing. Examples of applications using these buffers include storage and distribution of concentrated buffer, tangential flow filtration, pH re-equilibration, and final formulation.

The study will investigate the mixing performances for buffer preparation at different volumes of single-use bags (50 L, 200 L, 1,000 L), and at different transient concentrations of powders (16 g/L, 73 g/L, 224 g/L) before the final dilution.

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 following powders used in the buffer preparation:

- TRI buffer: Citric Acid, Sodium Citrate, Sodium Chloride (final concentration = 14.4 g/L)
- Tris buffer: Tris, Tris HCl, Sodium Chloride (final concentration = 67.6 g/L)
- 20x PBS (final concentration = 191 g/L)

The mixing times are determined by conductivity and visual inspection of the solution in the Flexel[®] Bag for Magnetic Mixer. The usual procedure for a buffer preparation includes before final dilution the incorporation of the powders in the bag partially filled with WFI. For our study, the bags were respectively filled to 80% of the nominal volume for the concentrated PBS and 90% for the TRI and Tris buffers.



200L Flexel[®] with Magnetic Mixer Technology

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

The actual concentrations of salts during the mixing phase before the final dilution are:

- TRI buffer: 16 g/L
(fluid expansion due to the powders is negligible)
- Tris buffer: 73 g/L (density = 1.046 kg/L)
- 20x PBS: 224 g/L (density = 1.162 kg/L)

Then the final step of the process consists in the addition of water to achieve the expected buffer concentrations.

Materials and methods

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

1. Standard Flexel® Bag for Magnetic Mixer (50L: FMB114867, 200L: FMB114893, 1,000L: FMB114896)
2. Powder transfer bag 15L (ref. FMA114008)
3. Pallettank for LevMixer® and Magnetic Mixer¹
(50 L: FXC110820, 200L: FXC110821, 1,000L: FXC113384)
4. Magnetic Mixer Drive Unit, 230V,
EU power cord (ref. LT-DU-006-EU)
5. Powder bag holder 200–400–650L (ref. FXA114344)
6. SAFC® powders with the following formulations:
TRI buffer (product number 44078 – dry powder packaging of 10 kg):
 - Citric acid anhydrous (0.04 g/L – 0.2 mM)
 - Sodium chloride (8.6 g/L – 147mM)
 - Sodium citrate, 2 H₂O (5.8 g/L – 20 mM)
 Tris buffer (product number 44077 – dry powder packaging of 10 kg):
 - Sodium chloride (58.44 g/L – 1M)
 - Tris (6.06 g/L – 50 mM)
 - Tris HCl (3.15 g/L – 20 mM)
 Concentrated 20x PBS (product number 44079 – dry powder packaging of 10 kg):
 - Potassium chloride (4 g/L – 54 mM)
 - Potassium phosphate, monobasic, anhydrous (4 g/L – 29 mM)
 - Sodium chloride (160 g/L – 2.7 M)
 - Sodium phosphate, dibasic, anhydrous (23 g/L – 162 mM).
7. Conductivity sensor: WTW InoLab Cond 740i

Method used:

1. The buffer is prepared in standard Flexel® Bags for Magnetic Mixer filled with deionised water to 80% of the final volume for the concentrated PBS, and 90% of the final volume for the Tris and TRI buffers.
2. Impeller speed is set to the maximum speed of 300 rpm to optimize the powders hydration and dispersion.
3. The powders are incorporated in the Flexel® Bags for Magnetic Mixer using either:
 - SAFC® packaging for dry powder,
 - or Sartorius Stedim Biotech 15 L powder transfer bag for a contained transfer to the mixing bag assembly.

4. Two mixing times are monitored from the addition of buffer powders:

- 4.1 "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 (see Fig. 1).

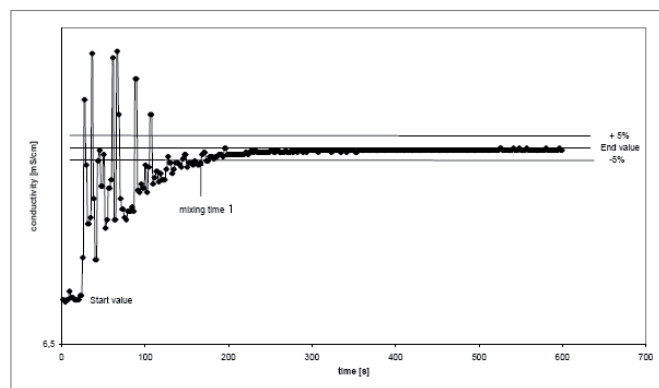


Fig. 1: General principle of mixing time determination via conductivity

- 4.2 "mixing time 2" is determined by a visual inspection.

The "mixing time 2" corresponds to the time when all suspended particles are visually dissolved.

Results and discussions

1. Mixing performances vs. volume of buffer preparation

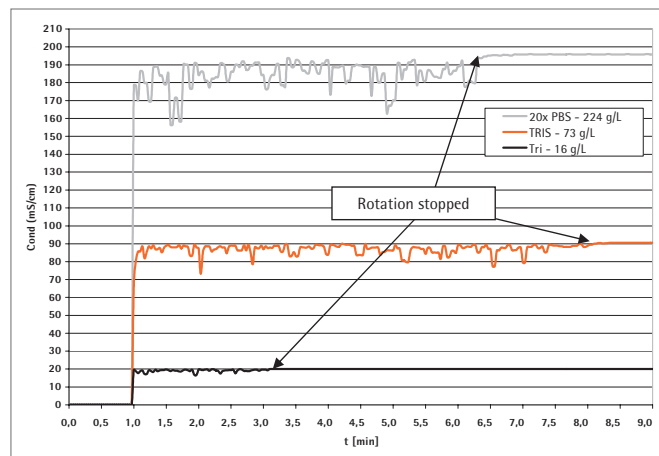


Fig. 2: Buffer preparations in 50L Flexel® Bag for Magnetic Mixer

50 L Flexel® Bag for Magnetic Mixer: the vortex at 300 rpm in the 50L 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. The mixing time is at least as good as with the larger scale.

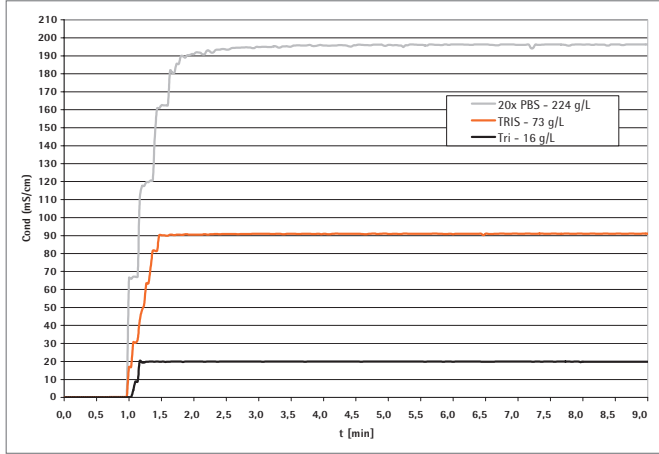


Fig. 3: Buffer preparations in 200L Flexel® Bag for Magnetic Mixer

200L Flexel® Bag for Magnetic Mixer: 4 SAFC® bucket liners were used to transfer 38.2 kg of PBS into the mixing bag as shown by the multiple steps on the conductivity signal for PBS buffer preparation (see Fig. 3). A single Sartorius Stedim Biotech powder transfer bag was used for the preparation of Tris (13.5 kg) and sodium citrate | citric acid (2.9 kg), which is reflected by continuous increase of conductivity signal. A stable plateau conductivity value was then quickly achieved for the three buffers, indicative of the mixing completion.

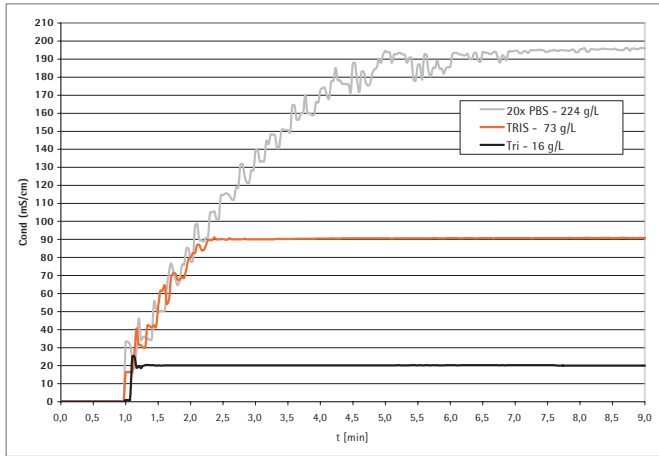


Fig. 4: Buffer preparations in 1,000L Flexel® Bag for Magnetic Mixer

1,000L Flexel® Bag for Magnetic Mixer: 19 SAFC® bucket liners were used to transfer 191 kg of PBS, 7 liners for the Tris (66.7 kg) and a single one for the sodium citrate | citric acid (14.4 kg). The multiple powder addition steps are visible on the conductivity graph (see Fig. 4). Again a stable value of conductivity was monitored at the mixing completion.

2. Mixing performances vs buffer type – determination of the mixing times

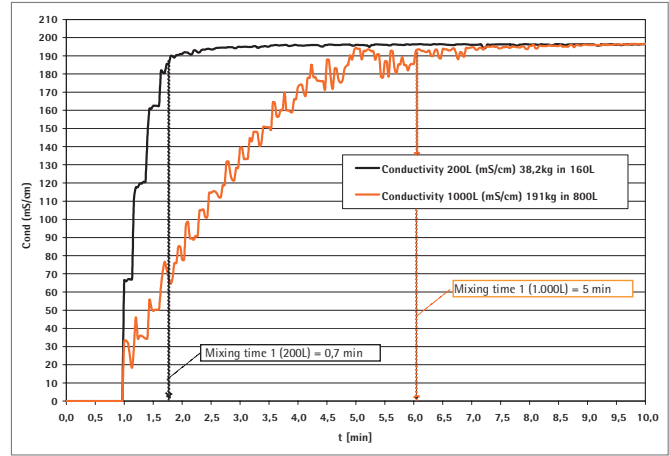


Fig. 5: PBS 20x buffer preparation in Flexel® Bag for Magnetic Mixer

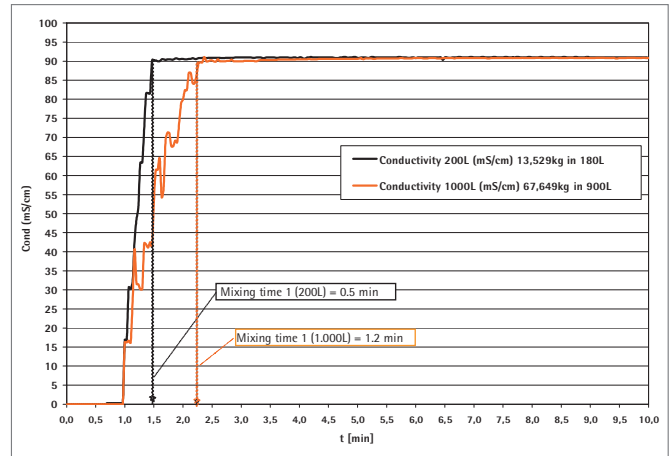


Fig. 6: TRIS buffer preparation in Flexel® Bag for Magnetic Mixer

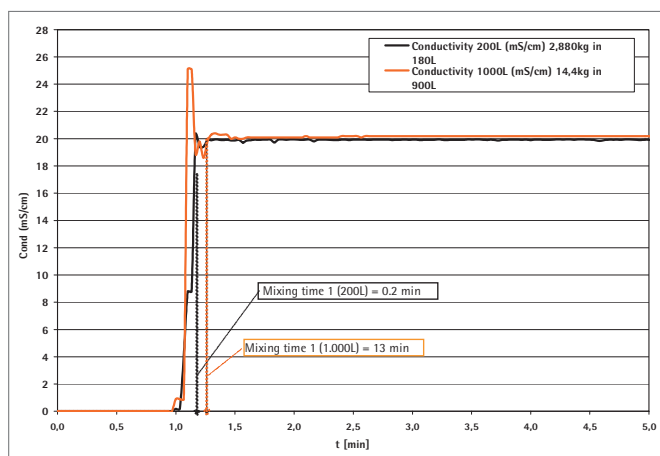


Fig. 7: Sodium citrate | citric acid/NaCl buffer preparation in Flexel® Bag for Magnetic Mixer

Bag Volume (L)		50	200	1000
Buffer powders				
Citric Acid Sodium Citrate NaCl (16.0 g/L)	Based on conductivity		0,2 min	0,3 min
	Based on visual inspection	<5 min	<5 min	<5 min
Tris Tris HCl NaCl (73.0 g/L)	Based on conductivity		0,5 min	1,2 min
	Based on visual inspection	<5 min	6 min	10 min
20x PBS (224.0 g/L)	Based on conductivity		0,7 min	5 min
	Based on visual inspection	13 min	14 min	20 min

Fig. 8: Overview on mixing times via conductivity and visual inspection

* conductivity values not stable at 50L due to presence of air bubbles around the conductivity cell (strong vortex)



200L Palletank for Magnetic Mixer equipped with the powder transfer bag for the mixing trial

General comments:

- The mixing times reported in this study include the transfer time of the multiple Sartorius Stedim Biotech powder transfer bags (for the 200 L scale experiment) or SAFC® bucket liners (50 and 1,000 L experiments) into the mixing bag assembly.
- A rapid dissolution of the powder salts was observed for the three buffers thanks to the strong mixing torque at 300 rpm (see Fig. 5-7).
- For the 200 L and 1,000 L volume, the conductivity of the solutions reaches a stable value in few minutes. However, some fine particulates of salt can still be visually observed in the solution. The agitation at 300 rpm was maintained until the solution became totally clear. This visual control is facilitated by the large windows of the Palletank.
- For all cases a 20 x increase in buffer preparation volumes (50L to 1,000 L) only leads to a 2x increase in the mixing times.
- The worst-case buffer preparation condition (1,000 L 20X PBS with mixing of 191 kg of salts in 800 L of water) resulted in a mixing time of 20 min.
- To the mixing times presented in Fig. 8, additional time for equipment set up and water filling to 80% as well as time for filling after mixing from 80 to 100% to reach final volume and concentration should be taken into consideration for total process time calculation.

Conclusion

- Large volume buffer solutions are quick and easy to prepare using the combination of ready to use buffer formulations and the high efficiency mixing of the Flexel® with Magnetic Mixer Technology.
- The contained processing conditions with the closed powder transfer bag docked onto the sterile Flexel® Bag for Magnetic Mixer are favourable to maintain low bioburden and to reduce the minimum exposure of the operator to chemicals.
- The platform provides a single-use scalable buffer preparation capability with a range of Flexel® Bags including volumes of 50L, 100 L, 200 L, 400 L, 650 L and 1,000 L.
- Flexel® for Magnetic Mixing system integrated in a FlexAct BP system provides monitoring and automation capability to better control all the operations of a buffer preparation step.

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
Modectcenterstrasse 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 Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
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