



sartorius stedim
biotech

Scalable buffer preparation with single-use LevMixer® technology



Application
Note

#14

#15

#16

#17

#18

turning science into solutions

Executive Summary

This application note presents the results of buffer preparations using the Flexel® for LevMixer®¹ technology. The method combines ready to use buffer formulations from SAFC®, with scalable, high efficiency, single use mixing systems provided by Sartorius Stedim Biotech.

The performances presented in this application note are achieved with the superconducting drive unit that enables a levitation of the impeller encapsulated into the Flexel® Bag. This technology is the alternative to the Flexel® for Magnetic Mixer² technology which is part 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 single-use mixing system enables a rapid dissolution and dispersion of the buffer powders in liquid for volumes of 50L to 1,000L. The performances of the single-use mixing system are characterized with quantitative (conductivity measurement) and qualitative (visual inspection) techniques.

As for the Flexel® for Magnetic Mixer technology, the LevMixer® technology provides seamless scale up and rapid mixing for process development and GMP manufacturing of buffers at 16g/L and 73 g/L. For the concentrated buffer at 224 g/L prepared in this study, the Flexel® for Magnetic Mixer technology is the preferred technology thanks to its higher rotation speed which minimizes the time for dissolution (refer to the related application note #14 for details).

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 used for this application study is the Flexel® for LevMixer® with volumes of 50L, 200L and 1000L. The levitated impeller enables a rotation speed up to 180 rpm, providing an efficient 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 at different volumes of single-use bags (50L, 200L, 1000L), and at different 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 LevMixer® technology to dissolve the following powders used in the buffers 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 LevMixer®.

The usual procedure for a buffer preparation requires the incorporation of the powders in the bag partially filled with WFI. For our study, the bags were respectively filled to 90% of the nominal bag's volume for the TRI and Tris buffer and 80% for the concentrated PBS.

The actual concentrations of salts for which mixing times are measured 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.

¹ LevMixer® is a trademark of Pall Corporation and this product uses Pall patented LevMixer® technology.

² This product uses Pall patented Magnetic Mixer technology.

All information on patents can be found at Pall.com/patents.



Materials and methods

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

1. Standard Flexel[®] Bag for LevMixer[®] (50L: FXB111567, 200L: FXB111420, 1000L: FXB111569)
2. Powder Transfer Bag 30L (ref. FMA114009)
3. Palletank for LevMixer[®] and Magnetic Mixer¹ (50L: FXC110820, 200L: FXC110821, 1000L: FXC113384)
4. Superconducting drive unit, 230V, EU power cord (ref. LT-DBTL-007)
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 10kg):
 - Citric acid anhydrous (0.04 g/L – 0.2mM)
 - Sodium chloride (8.6 g/L – 147mM)
 - Sodium citrate, 2 H₂O (5.8 g/L – 20mM)
 Tris buffer (product number 44077 – dry powder packaging of 10kg):
 - Sodium chloride (58.44 g/L – 1M)
 - Tris (6.06 g/L – 50mM)
 - Tris HCl (3.15 g/L – 20mM)
 Concentrated 20x PBS (product number 44079 – dry powder packaging of 10kg):
 - Potassium chloride (4 g/L – 54mM)
 - Potassium phosphate, monobasic, anhydrous (4 g/L – 29mM)
 - Sodium chloride (160 g/L – 2.7M)
 - Sodium phosphate, dibasic, anhydrous (23 g/L – 162mM).
7. Conductivity sensor: WTW InoLab Cond 740i

Method used:

1. The buffer is prepared in standard Flexel[®] Bag for LevMixer[®] 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 180 rpm to optimize the powders hydration and dispersion.
3. The powders are incorporated in the Flexel[®] Bag for LevMixer[®] using either:
 - SAFC[®] packaging for dry powder,
 - or Sartorius Stedim Biotech 30L 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.

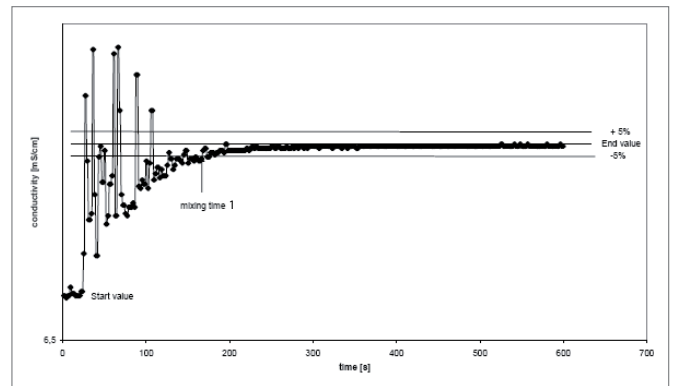


Fig. 1: General principle of mixing time determination via conductivity

- 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

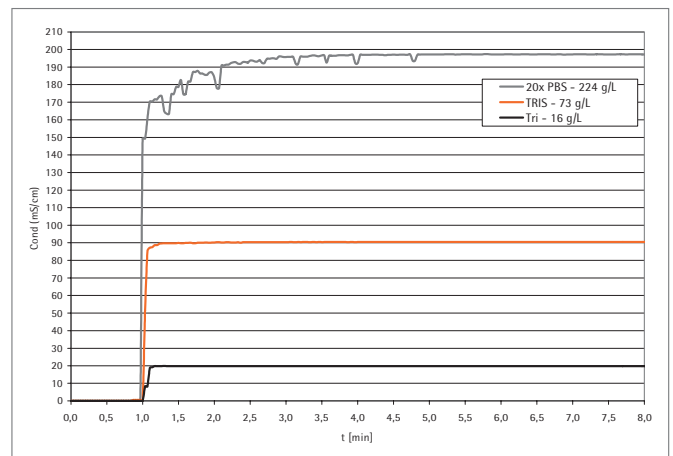


Fig. 2: Buffer preparations in 50L Flexel[®] Bag for LevMixer[®]

50L Flexel[®] Bag for LevMixer[®]: A single bucket liner from SAFC[®] was used to transfer the powders for each of the 3 buffers, thus explaining the continuous conductivity increase. A stable value was then quickly observed, indicative of the mixing completion.

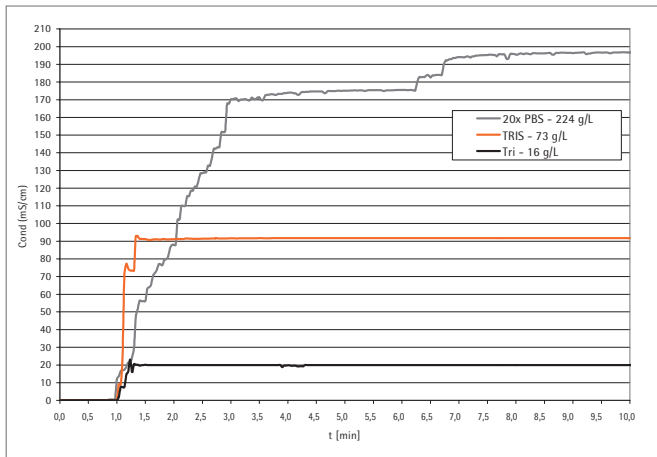


Fig. 3: Buffer preparations in 200L Flexel® Bag for LevMixer®

200L Flexel® Bag for LevMixer®: 2 separate Sartorius Stedim Bio-
tech 30L powder transfer bags were used to add the 38.3 kilos of
PBS. The time for disconnection and connection of the 2nd
powder transfer bag explains the stable period of conductivity
observed between the 3rd and the 6th minute during the PBS
addition. Two SAFC® bucket liners were used for the preparation
of TRIS (13.5 kg) and a single one for sodium citrate|citric acid
(2.9 kg), which is reflected by continuous increase of conductiv-
ity signal. A stable plateau conductivity value was then quickly
achieved for the three buffers, indicative of the mixing completion.

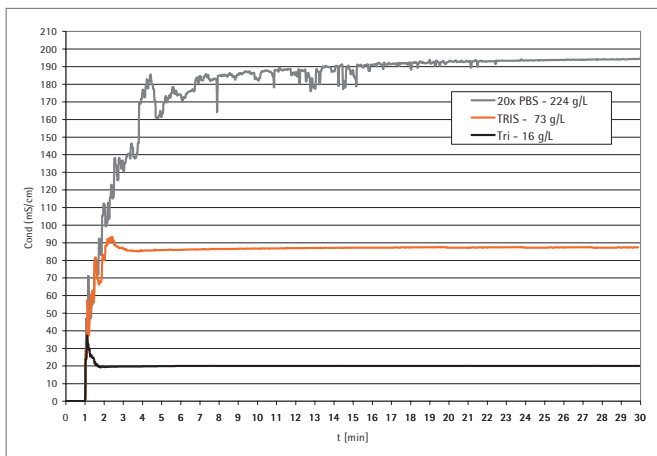


Fig. 4: Buffer preparations in 1000L Flexel® Bag for LevMixer®

1000L Flexel® Bag for LevMixer®: 19 SAFC® bucket liners were
used to transfer 191kg of PBS, 7 liners for the Tris (66.7kg) and a
single one for the sodium citrate|citric acid (14.4kg). The multiple
powder addition steps are visible on the conductivity graph.
Again a stable value of conductivity was monitored at the mixing
completion.

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

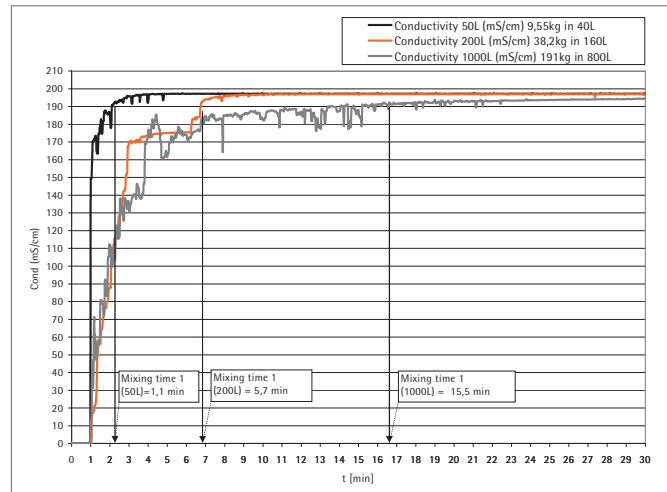


Fig. 5: PBS 20x buffer preparation in Flexel® Bag for LevMixer®

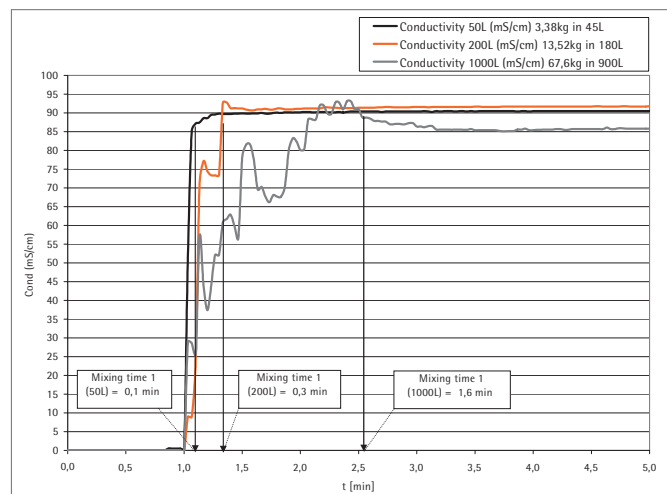


Fig. 6: TRIS buffer preparation in Flexel® Bag for LevMixer®

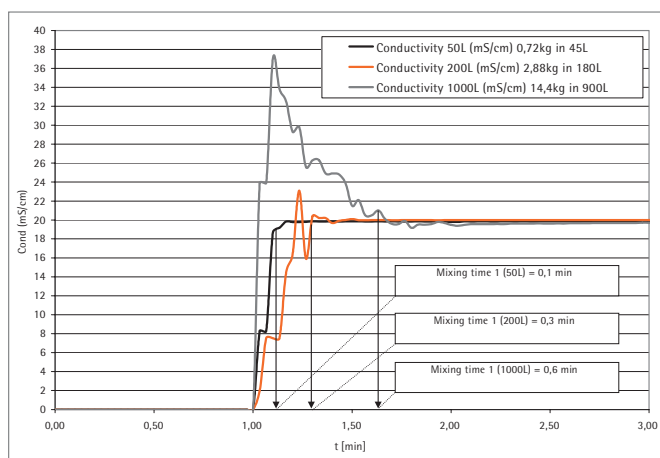


Fig. 7: Sodium citrate | citric acid | NaCl buffer preparation in Flexel® Bag for LevMixer®

Bag Volume (L)		50	200	1000
Powders Nature				
Citric Acid Sodium Citrate NaCl (16.0 g/L)	Based on conductivity	< 1 min	< 1 min	< 1 min
	Based on visual inspection	< 5 min	< 5 min	< 5 min
Tris Tris HCl NaCl (73.0 g/L)	Based on conductivity	< 1 min	< 1 min	1-2 min
	Based on visual inspection	< 5 min	8 min	14 min
20x PBS (224.0 g/L)	Based on conductivity	1-2 min	5-6 min ^(*)	15-16 min
	Based on visual inspection	10 min	32 min	74 min

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

* 5-6 minutes including 3 minutes for Sartorius Stedim Biotech's exchange of powder bag.

General comments:

- The mixing times reported in this study include the transfer time of the multiple Sartorius Stedim Biotech Powder Transfer Bag (for the 200L scale experiment) or SAFC® bucket liners (50, 200 and 1000L experiments) into the mixing bag assembly.
- For each volume of Flexel® Bag, the conductivity of the solutions reaches a stable value in a few minutes. However, some fine particles of salt can still be visually observed in the solution. The agitation at 180 rpm was maintained until the solution became totally clear. This visual control is facilitated by the large windows of the Palletank.
- With respective mixing times below 2 minutes (by conductivity) or 14 minutes (by visual observation), a rapid dissolution of the salts was observed with:
 - the TRIS (50L, 200L, 1000L)
 - the Sodium Citrate, Citric Acid (50L, 200L, 1000L)
 - the 20x PBS (50L).
- The worst-case conditions identified with the 20x PBS in the 200L and 1000L Flexel® Bags for LevMixer® require extended times for complete powder dissolutions. Nevertheless a stable value of the conductivity and a complete dissolution were achieved like with the lower concentrations and volumes.

To the mixing time 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 efficient mixing of the Flexel® with LevMixer® Technology. For the concentrated buffer 20x PBS prepared in large volumes (200L and 1000L), the Magnetic Mixer Technology is recommended to optimize the mixing times thanks to the higher rotation speed (300 rpm with Magnetic Mixer versus 180 rpm with LevMixer®).
- The contained processing conditions with the closed Powder Transfer Bag docked onto the sterile Flexel® Bag for LevMixer® 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 buffer preparation capability with a range of Flexel® Bags including volumes of 50L, 100L, 200L, 400L, 650L and 1000L.

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