

# 1X PBS (Phosphate Buffered Saline) Buffer Mixing Optimization

## Using the Mobius® Single-Use Mixing Solution with Process Chemicals and Filters

### Introduction

This study develops optimal process conditions for mixing and filtering of 1X PBS buffer using the 50 L, 200 L, and 1000 L Mobius® Single-Use Mixing Solution, Durapore® and Millipore Express® filters.

The Mobius® Single-use Mixing Solution delivers advanced technology for mixing pharmaceutical ingredients from intermediate to final drug products and for the preparation of process solutions, such as buffers and media. Each single-use mixing container includes a magnetically-driven levitating impeller for improved mixing consistency and efficiency. The Mobius® Single-use Mixing Solution includes a powder delivery system, 10 L, 50 L, 100 L, 200 L, 500 L, and 1000 L single-use mixers and the Mobius® FlexReady Solution for Buffer and Media Preparation.

Durapore® 0.22 µm, Millipore Express® SHC (0.5/0.2 µm), and Millipore Express® SHF (0.2 µm) sterilizing-grade filters were used in this study. Durapore® filters contain polyvinylidene fluoride (PVDF) membrane and Millipore Express® filters contain polyethersulfone (PES) membrane. These sterilizing-grade membranes are offered in capsule and cartridge formats in a variety of pore sizes to meet different sterility assurance needs. The capsules are gamma and autoclave compatible and can be purchased gamma sterilized. The gamma compatible capsules are incorporated directly into the disposable Flexware® assemblies of the Mobius® FlexReady Solution for Buffer and Media Preparation, resulting in a sterile, fully-closed flowpath.

Many of our buffer products bear the Emprove® Expert and Emprove® Essential trademark. The Emprove® product lines have been developed to make work safer, faster and more cost-efficient, and offer regulatory documentation that strengthens the quality section of registration dossiers.

### Experimental Methods

#### Mixing Study

Three batches of 1X PBS were prepared at 50 L, 200 L, and 1000 L volumes using the following methodology:

1. The bag was installed into the mixer container and inflated prior to use.
2. The scale/load cells were tared with the empty mixer vessel. A floor scale was used for the 50 L and 200 L mixer vessels. The 1,000 L vessel comes with on-board load cells.
3. Water was added to approximately 90% of the final target volume.
4. The mixer vessel was turned on and the impeller was set to the maximum setting:
  - Mobius® MIX50: 1000 RPM
  - Mobius® MIX200: 500 RPM
  - Mobius® MIX1000: 250 RPM
5. The Emprove® Expert chemicals used to prepare 1X PBS buffer were weighed into a powder delivery bag, the bag was connected to the mixer bag and the powder delivery valve was opened to transfer the powder to the mixer vessel as one bulk addition. For the 1,000 L scale preparation, the Mobius® hoist was used to lift the powder bag above the mix vessel and hold it in place while the powder bag was connected to the mix bag. The hoist was not needed to connect the powder bag to the 50 L and 200 L mixer bags. **Table 1** details the chemical materials needed to prepare individual batches of PBS, including the final concentration and final total powder masses.

- Samples were taken from the top of the mixer vessel and from the needle-free sample ports beginning immediately after adding the powder into the mixer vessel. Samples were returned back to the mixer vessel after each measurement.
- The pH and conductivity of the buffer was measured every 2 minutes. When the measurements at the top and bottom were equal and stable for 10 minutes, the buffer was considered mixed. At this point water was added to 100% of the final target with the impeller still running.

**Table 1. Emprove® Expert Chemical Requirements for Individual Batches of 1X PBS buffer**

Chemical	Conc [g/L]	Weight [g]		
		50 L	200 L	1000 L
Sodium Chloride	8	400	1600	8000
di-Sodium Hydrogenphosphate, Dihydrate	1.78	89	356	1780
Potassium Dihydrogen Phosphate	0.27	14	54	270
<b>Total Powder Weight [kg]</b>	<b>N/A</b>	<b>0.503</b>	<b>2.01</b>	<b>10.05</b>

## Filtration Study

After the 1X PBS buffer mixing study was performed at 50 L scale, the filterability of the buffer was tested using the Vmax™ method at a constant pressure of 10 psi to determine which filter type and area would be most appropriate for the three different batch sizes. The Vmax™ method is an accelerated screening test based on the gradual pore size plugging model, and estimates scale up filter area requirements. Three replicates were performed for each filter to ensure data was repeatable and reliable. OptiScale® 25 devices (3.5 cm<sup>2</sup> membrane area) were used during Vmax™ testing. **Table 2** outlines the filters that were evaluated in this study.

**Table 2. Filter Evaluation**

Filter	Chemistry	Pore Size
Millipore Express® SHC	Hydrophilic PES	0.5/ 0.2 µm
Millipore Express® SHF	Hydrophilic PES	0.2 µm
Durapore®	Hydrophilic PVDF	0.22 µm

## Results and Discussion

### Mixing Study

**Table 3** details the mixing times required to get the first stable pH and conductivity readings prior to filling to 100% of the final volume (i.e. when the readings of samples taken from the top of the carrier and from the sample ports matched each other). They were 12, 6 and 10 minutes for the 50 L, 200 L, and 1000 L scales, respectively. This demonstrates that the mixer technology is effective and efficient across a wide range of volumes. Once the final water addition is complete, it is recommended to continue mixing and verify that the pH and conductivity values remain stable for a minimum of 10 minutes before filtering or using the buffer.

**Table 3. Mixing Time Requirements for 1X PBS Buffer Preparation at 50 L, 200 L, and 1000 L Scales**

Step	Step Duration		
	50 L	200 L	1000 L
PBS buffer addition	< 3 seconds	< 3 seconds	9 seconds
Mixing time to get the first stable value of pH and conductivity	12 minutes	6 minutes	10 minutes

In this study no attempt was made to match the mixing times of the two different scales of preparation by adjusting the impeller speed or initial fill volume. Instead, the maximum impeller speed available at each scale was used with a fixed initial volume based on experience and ease-of-use as described above. Separate studies have been performed with the Mobius® single-use mixing products to determine a scaling algorithm for both floating and sinking powders. Please contact your local representative for more information.

## Filtration Study

The optimal filter choice and area requirement for sterile filtration of the 1X PBS buffer was determined using the Vmax™ method at a constant pressure of 10 psi. This method is an accelerated screening test based on the gradual pore size plugging model and it estimates scale up filter area requirements. Vmax™ testing was performed in triplicate for all filters shown in **table 2** using the freshly prepared 50 L batch of 1X PBS buffer. Since PBS buffer is non-plugging, the Vmax™ results were then used to project filtration area requirements for a range of batch sizes and process times using the following equation:

$$\text{Equation 1: } A_{\text{min}} = V_B / (J_{\text{mean}} * t_B)$$

Where:

A<sub>min</sub>: minimum membrane area of the filter [m<sup>2</sup>]  
V<sub>B</sub>: batch volume [L]  
J<sub>mean</sub>: mean flux [L/ (m<sup>2</sup>.h)]  
t<sub>B</sub>: batch process time [h]

In addition, the K<sub>Buf</sub> method of determining sterile filter requirements for non-plugging buffers was also used and these results were compared to the filter recommendations generated by the Vmax™ method. K<sub>Buf</sub> is an efficient and simplified filter sizing tool for buffer filtration which is based on well characterized membrane permeability values. The K<sub>Buf</sub> sizing method has been shown to closely approximate sterile filter sizing requirements for buffer filtration applications generated through the more laborious Vmax™ sizing method. Filter sizing via the K<sub>Buf</sub> method is determined using the following equation:

$$\# \text{ of } 10'' \text{ Filter Cartridges} = V_B \times K_{\text{Buf}} / t_B \times \Delta P$$

Where:

V<sub>B</sub>: batch volume [L]  
K<sub>Buf</sub>: sizing constant for a specific filter type which can be applied to water-like (viscosity = 1cP), non-plugging buffers  
t<sub>B</sub>: batch filtration time [h]  
ΔP: filtration process differential pressure [psid]

In all cases, the filter sizing recommendations resulting from the Vmax™ and the K<sub>Buf</sub> methods were within 10% of each other, supporting the continued use of K<sub>Buf</sub> as a simpler alternative sizing method for non-plugging buffers.

Of the three 0.2 μm filters evaluated in this study, Millipore Express® SHF membrane exhibited the best overall performance, requiring the lowest filter area. Millipore Express® SHC membrane required approximately 33% additional filter area and Durapore® 0.22 μm membrane required approximately 155% additional area to process the same batch volume in equivalent time.

**Table 4** summarizes the optimal filter capsule recommendations for 0.2 μm sterile filtration of 1X PBS buffer batches ranging from 10 L to 1,000 L. Since the sizing calculation includes batch time, times were selected that seemed reasonable and would be achievable using the pumping system on the Mobius® FlexReady Solution for Buffer and Media Filtration, which can operate at flow rates up to 10 L/min. In all cases, the final filter products chosen are gamma-compatible capsules that can be incorporated directly into Flexware® single-use assemblies, resulting in sterile, fully-closed flowpaths for the final buffer processing. In keeping with best practice guidance, a minimum 50% safety factor was applied to all minimum areas and the filter capsules with the closest membrane area at or above this safety factor were selected.

**Table 4. Recommendations for sterile filtration of 10–1,000 L batch sizes of 1X PBS buffer**

Batch Size [L]	Filtration Time [min]	Recommended Filter
10	10	Opticap® XL 150 with Millipore Express® SHF membrane
50	15	Opticap® XL 300 with Millipore Express® SHF membrane
100	20	Opticap® XL 600 with Millipore Express® SHF membrane
200	30	Opticap® XL 600 with Millipore Express® SHF membrane
500	60	Opticap® XL 600 with Millipore Express® SHF membrane
1000	120	Opticap® XL 600 with Millipore Express® SHF membrane

Since 1X PBS buffer is a non-plugging solution, filter sizing is dependent on flux and batch time, not volumetric capacity. Therefore, the selected filtration time determines the required membrane area. If a smaller device is desired, this can be achieved by increasing the batch filtration time.

Ultimately, using the recommendations shown above, the Mobius® FlexReady Solution for Buffer and Media Filtration was successfully used to sterile filter both the 200 L and 1,000 L batches of 1X PBS buffer that were prepared in the Mobius® Single-use Mixers.

## Conclusions

This study highlights an efficient way to prepare and filter 50 L, 200 L, and 1000 L batches of 1X PBS buffer using Mobius® FlexReady Solution for Buffer and Media Preparation, Mobius® Single-Use Mixers and sterile filters. Millipore Express® SHF filters were found to be the optimal choice for sterile filtration of the 1X PBS buffer for all batch sizes.

### Best Practices for Mixing PBS

- It is recommended to inflate the mixer bags with air prior to the addition of water to ensure proper alignment with the mix vessel.
- The clamp on the bottom outlet port should be placed as close as possible to the bag to minimize dead leg volume where salts could accumulate.
- Filling the mixer with water to approximately 90% of the target final volume prior to adding powders is recommended.
- In order to achieve lowest mix times, it is recommended that the impeller be set at the maximum RPM setting. However, if scalable mix times is a concern, contact your local representative for information.
- The chemical components should be added to the mix bag at a rate which prevents the accumulation of large amounts of powder at the bottom of the mix bag, in order to avoid settling in the outlet port.
- Once all of the components are visually dissolved, the pH can be adjusted, if necessary. Finally the solution can be brought to 100% of final target volume by adding water with the impeller still running.

MilliporeSigma  
400 Summit Drive  
Burlington, MA 01803

## To place an order or receive technical assistance

In the U.S. and Canada, call toll-free 1-800-645-5476

For other countries across Europe and the world, please visit: [EMDMillipore.com/offices](https://www.emdmillipore.com/offices)

For Technical Service, please visit: [EMDMillipore.com/techservice](https://www.emdmillipore.com/techservice)

[EMDMillipore.com](https://www.emdmillipore.com)

