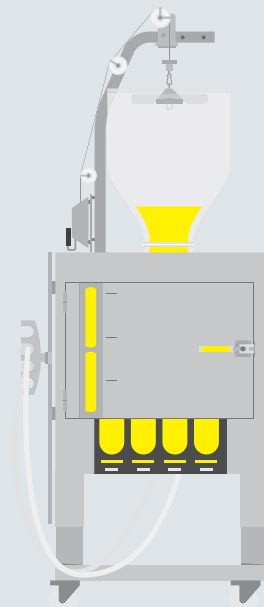


Ready-to-Hydrate Buffers and Media

Pre-Weighed Powders in Flexel® Tank Liners and Powder Transfer Bags for Media and Buffer Preparation Applications



Benefits

Reducing Time and Contamination Risks

Powder Bags are connectable to any mixing system via their tri-clamp connector.

Powders are manufactured according to the highest quality standards from qualified, compendial grade raw materials.

Available for Sartorius proprietary media as well as any custom process formulations.

Consistency From Batch to Batch

Powders are manufactured in ISO8 clean-rooms following cGMP referentials according to users' unique specifications.

Ready-to-Hydrate Buffers and Media

Once hydrated, only sterile filtration is left to do. If required, pH adjustment can be easily achieved with our off-the-shelf sterile NaOH 1M in Flexsafe® bag.

Product Information

Powder Buffers and Media Prefilled Bags are pre-weighed and designed to match requirements for the smaller volumes up to manufacturing scale. Flexel® Tank Liners and Powder bags will respectively match open and closed mixing solutions.

As media and buffers composition vary from one process to another, our team of experts will provide support for the evaluation of any custom request.

Applications

Our manufacturing capabilities cover all powder compositions from single component products to complex blends; ranging from upstream cell culture media to downstream purification process formulations. Two primary packaging options are available in order to match with any mixing systems.

Powder Prefilled Bags are single-use primary packaging solutions which reduce risks of contamination by avoiding dust when preparing large powder volumes.

In bioprocessing, **the safety of the operators** is a critical aspect increasingly outlined by regulatory agencies. This is where **Ready-to-Hydrate Powders help improve environment conditions** (including ATEX) as they offer a dust-free contained powder handling for mixing.

From regular buckets to drums, to pre-filled bags, get each of your powder formulations in your preferred primary packaging.

Technical Data

Specifications

Features	Benefits	Comment
Dust-free	Reduce risks of contamination	4" or 8" Tri-Clamp connection for Powder Transfer Bags
Non-Animal Origin	Regulatory friendly for research or further manufacturing use only	Certificate available on request*
Precise Filling	One bag, one prep. Reducing errors and batch to batch variability	Weighing tolerance ±1%
Milled powder	Easier and faster powder dissolution	Homogeneous particle size distribution
Robust packaging	Suitable for GMP operations	Each Prefilled bag is double-wrapped in vacuum-sealed pouches

* Powder Buffers and Chemically Defined Media are non-animal origin certified. Powders are manufactured according to cGMP referentials in ISO-certified facilities.

Characteristics

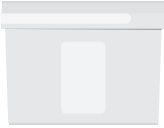



Sartorius' range of WFI-Quality Water in Flexsafe® bags are recommended for the powder hydration step which can be achieved quickly and easily even in worse-case conditions with Flexsafe® Pro Mixer high mixing performances.

Ready-to-Hydrate Powders require minimal setup and reduce footprint for the media and buffer preparation. The hydration process is achieved in a few minutes only; the milling process for the powder manufacturing ensures a homogeneous particle size distribution.

The joined knowledge and expertise in powder and film manufacturing ensures the supply of a reliable and convenient solution.

Sartopore® 2 filters are recommended for the sterilization-grade filtration step.

Complete Range of Packaging Solutions for Powder Media and Buffers

		10 g to 2 kg	5 kg to 20 kg	500 g to 10kg	2 kg to 20 kg	
Bucket		■				
Drum			■			
Flexel® Tankliner				■		Suitable for Open Mixing Systems
Powder Transfer Bag					■	Suitable for closed mixing systems including Flexsafe® ProMixer
				Up to 200 L	Up to 2,000 L*	

* Based on a regular PBS formulation ~10 g/L

Ordering Information*

Powder Buffers Prefilled Bags are also available as custom products. They will be manufactured to the exact specifications requested and delivered with the appropriate documentation. Contact your local Account Manager or Sartorius expert for more information. Below are product information for our standard, off-the-shelf powder Phosphate Buffer Saline (PBS) references.

Product Description	Product	Packaging	Volume	Order Number
Ready-to-Hydrate Buffer	PBS	Flexel® Liner	50 L	CQB3FA0001
Ready-to-Hydrate Buffer	PBS	Flexel® Liner	100 L	CQB3FA0002
Ready-to-Hydrate Buffer	PBS	Flexel® Liner	200 L	CQB3FA0003
Ready-to-Hydrate Buffer	PBS	Powder Transfer Bag	100 L - in 15 L bag size	CQB3FA0004
Ready-to-Hydrate Buffer	PBS	Powder Transfer Bag	200 L - in 15 L bag size	CQB3FA0005
4"- 8" Tri-Clamp Connector				FMA114007

Additional Product Information

- Shelf life: 2 years
- Shipping and storage temperature: 2 °C to 30 °C
- Documentation:
Certificate of Analysis, Safety Datasheet

Standard PBS pH 7.4

Component	g/L	CAS #	Grade
KH ₂ PO ₄	0.283	7778-77-0	EP USP
Na ₂ HPO ₄	1.124	7558-79-4	EP USP
NaCl	9.000	7647-14-5	EP USP

Complementary Products

Product Description	Product	Packaging	Volume	Order Number
WFI-Quality Water	WFI	Flexsafe®	20 L (EMEA and Asia)	CFB3FA3006
WFI-Quality Water	WFI	Flexsafe®	20 L (Americas)	CFB3FA3106
WFI-Quality Water	WFI	Flexsafe®	200 L (EMEA and Asia)	CFB3FA3009
WFI-Quality Water	WFI	Flexsafe®	20 L (Americas)	CFB3FA3109
NaOH 1M	Sodium Hydroxide	Flexsafe®	20 L	CFB3FA7006


* Other product formats and formulations available. For more information please contact: Europe | Asia: CellCultureMedia.EU@sartorius.com, Americas: CellCultureMedia.NA@sartorius.com

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com