

Sodium Hydroxide

NaOH 1M in
20 L Flexsafe® Bag



Features & Benefits

- Highly concentrated solutions can help reduce footprint and warehousing when diluted at the point of use
- Proven chemical compatibility with Flexsafe® film
- One single vendor to qualify for the sourcing of buffers & dangerous goods

Robust supply chain	Compendial grade raw materials
Regulatory-friendly for bioprocessing	Produced according to cGMP referentials
Sterile-grade filtration before filling	Sartopore® 2 filter family
Consistent & qualified process	WFI-quality water (EP) for liquid manufacturing
Fully documented	<ul style="list-style-type: none"> ▪ MSDS for first shipment or on request ▪ CoA provided with each batch

Highest Quality Standards

Content of the certificate of conformity

Test Method	QC-Release Parameter Specification
Appearance	Clear, colorless liquid
Density at 20 °C [g/cm ³]	For information
Conductivity [mS/cm]	For information
Assay [mol/L]	0.95 - 1.05
Test of sterility	Confirms
Shelf life	24 months

One single source for all bioprocessing buffers ranging from regular downstream buffers to Acids, Bases, Alcohols & Detergents

Applications

- Dangerous formulations often are required for cleaning, sanitization and specific downstream steps such as chromatography and virus inactivation

NaOH 1M is used for a broad variety of applications in the bioprocessing environment for 1M or lower concentrations. 20L Flexsafe® bags are the most appropriate format and a cGMP manufacturing process using WFI-quality water ensure the highest level of quality for a product that can be used for the following applications:

- Chromatography
 - flushing & activation of the membranes
 - packing and cleaning of the protein A
- Cleaning of reusable | stainless steel equipment
- Virus inactivation

Dedicated Packaging

Formulations used in downstream applications as well as for some specific process steps may be classified as dangerous goods; consequently dedicated labeling and packaging should be applied.

	10L - 20 L	50 L - 100 L - 200 L
Flexsafe® 2D bags	■	
Flexsafe® 3D bags for drums		■
UN-certified cardboard boxes	■	qualified for ADR Hazmat shipment worldwide
UN-certified drums		■ qualified for ADR Hazmat shipment worldwide

Each Flexsafe® 2D bag is wrapped in a sealed EAE* pouch.

* EAE is a coextruded multilayer film with a structure PE/PA/PE suitable for packaging medical devices

Ordering Information

Product Description	Packaging	Volume	Order Number
NaOH 1M	Flexsafe®	20 L	CFB3FA7006

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