

Thermo Scientific Nalgene Syringe Filters with Nylon Membrane

Cat. No.	Syringe Filter Description	Pore size, μm	Material	Packaging	No. per pack/ No. per case
195-2520 [†]	Sterile filter	0.2	Polypropylene	Ind. blister pack	50
195-2545 [†]	Sterile filter	0.45	Polypropylene	Ind. blister pack	50
196-2020	Non-sterile filter	0.2	Polypropylene	Bulk pack	50/300
196-2045	Non-sterile filter	0.45	Polypropylene	Bulk pack	50/300

[†]Certified non-pyrogenic and non-cytotoxic.

N/A – not applicable

These Nalgene[®] syringe filters are resistant to a wide range of solvents. Ideal for filtering samples and mobile phases used in gas chromatography and high pressure liquid chromatography. The Luer-Lok[®] design accommodates both standard and threaded Luer-Lok syringes. Nylon membrane is inherently

hydrophilic, contains no plasticizers or wetting agents and has low levels of extractables. Units are especially useful in critical applications requiring low contamination levels. Excellent for use with DMSO. Autoclavable.

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

Instructions for use

For either filtration method:

- For sterile filtration, always use a sterile syringe
- Use aseptic technique when required
- Always dispose of syringe and needle properly

Pressure Filtration

1. Fill syringe with sample (see "Caution," below)
2. Carefully remove the syringe filter from its package. Secure the syringe to the filter with a clockwise twist.
3. Hold the pre-filled syringe and filter vertically. This allows proper wetting of the membrane and helps prevent the formation of air blocks.
4. To begin filtration, press the syringe plunger.
5. The syringe filter may be used until resistance to pressure indicates clogging of the filter.

Caution: Use caution with syringes smaller than 10cc (mL). The pressure generated in these syringes may exceed the pressure limit of the syringe filter.

Aspiration or Vacuum Filtration

1. Carefully remove the syringe filter from its package. Secure the syringe to the filter with a clockwise twist. Attach the hub of the syringe needle to the outlet side of the syringe filter.
2. To filter liquid, retract the syringe plunger.
3. After filtration, remove the needle and syringe filter. This prevents particles trapped on the membrane from recontaminating the filtrate. Dispense filtrate by pressing filter syringe.

Nalgene Filterware is for research use only. It is not for *in vitro* diagnosis or parenterals.

WARNING:

Misuse of Nalgene products can be potentially dangerous. Before using this product, please refer to the appropriate Nalgene catalogs/inserts and the various warnings, information, instructions and chemical resistance charts. If any doubt exists about a specific use of Nalgene products, please contact Technical Service at 1-585-586-8800 or Fax to: 1-800-625-4363.
E-mail at: Technical.nalgene@thermofisher.com.

Pore size (µm)	Max. oper. pressure (psig/bar)	Surface area, cm ²	Bubble point (psig/bar) after air purge	Hold-up volume, mL	Nom. flow rate for water at (mL/min @ 14.7 psig=1 bar)
0.2	90/6.2	2.8	42/2.8	<0.15	35
0.45	90/6.2	2.8	25/1.7	<0.15	95

Chemical Resistance

These syringe filters are resistant to alcohols, aldehydes, ketones, esters, formamides, acetamides, dilute solutions of phenol or formaldehyde, strong alkalies, chlorinated or aromatic hydrocarbons, amines, freons, dilute hydrogen peroxide and organic acids.

Do not use with strong inorganic acids, such as concentrated hydrochloric acid, nitric, sulfuric or phosphoric acid; formic acid over 40%; acetic acid over 80%; concentrated phenols or cresols.

Autoclaving Guidelines

To autoclave, wrap syringe filter in kraft paper. Autoclave at 121°C/15 psig (1.03 bar) for 15-20 minutes. Store in a clean, dry place.

©2010 Thermo Fisher Scientific Inc. All rights reserved. Luer-Lok is a registered Trademark of Becton-Dickinson Co. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.thermoscientific.com

Asia: China Toll-free: 800-810-5118 or 400-650-5118; India: +91 22 6716 2200, India Toll-free: 1 800 22 8374;

Japan: +81 3 3816 3355 Other Asian countries: 65 68729717

Europe: Austria: +43 1 801 40 0; Belgium: +32 53 73 42 41; Denmark: +45 4631 2000; France: +33 2 2803 2180; Germany: +49 6184 90 6940, Germany Toll-free: 08001-536 376; Italy: +39 02 02 95059 or 434-254-375;

Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 100;

Russia/CIS: +7 (812) 703 42 15; Spain/Portugal: +34 93 223 09 18; Switzerland: +41 44 454 12 12;

UK/Ireland: +44 870 609 9203

North America: USA/Canada +1 585 586 8800; USA Toll-free: 800 625 4327

South America: USA sales support: +1 585 899 7198

Countries not listed: +49 6184 90 6940 or +33 2 2803 2180

8-0402-35 0310

Thermo
SCIENTIFIC