

Thermo Scientific Nalgene Syringe Filters with PES Membrane

**Cat. No. 180-1320 13 mm dia., 0.2 µm pore size,
acrylic housing**

**Cat. No. 194-2520 25 mm dia., 0.2 µm pore size,
acrylic housing**

**Cat. No. 194-2545 25 mm dia., 0.45 µm pore size,
acrylic housing**

Packaged individually, 50 per case

Nalgene® syringe filters with PES (polyethersulfone) membrane are specially designed to sterilize small volumes of aqueous liquids using a syringe. PES membranes offer lower protein binding, lower extractables, no wetting agents and higher flow rates than cellulosic membranes. They can be used with both standard slip-fit and Luer-Lok® syringes and accept standard hypodermic needles.

Nalgene PES membranes and syringe filters have good chemical resistance to aqueous solutions, glycols and alcohols. They are not recommended for use with esters, ethers, ketones, strong acids, or aliphatic, aromatic and halogenated hydrocarbons or their derivatives.

Nalgene syringe filters are preassembled, disposable and certified sterile, non-cytotoxic and non-pyrogenic.

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

Thermo Scientific Nalgene Syringe Filters with PES Membrane

Instructions for Use

Sterile Filtration

1. Always use a sterile syringe.
2. Use aseptic technique when required.
3. 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 inlet side of the filter with a clockwise twist.
3. Hold the pre-filled syringe and filter vertically. This allows proper wetting of the membrane and helps to prevent formation of air blocks.
4. Press syringe plunger to begin filtration.
5. The syringe filter may be used until resistance to pressure indicates clogging of the filter.

CAUTION: Use caution with syringes smaller than 10cc (10ml). The pressure generated in these syringes may exceed the 75 psig limit of the syringe filters.

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 syringe plunger.

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 any Thermo Scientific Nalgene product, please contact Technical Support at technical.nalgene@thermofisher.com or call 800.652.4327.

Specifications

Thermo Scientific Nalgene Syringe Filters with PES Membranes

Pore Size, μm	Max. Oper. Pressure (psig/bar)	Surface area, cm^2 (membrane dia.)		Hold-up volume (membrane dia.)	
		13mm	25mm	13mm	25mm
0.2	75/5.1	0.8	5.3	20 μl	0.15ml
0.45	75/5.1	NA	5.3	NA	0.15ml

Pore size, μm	Bubble Point (psig/bar)	Nom. Flow Rate (ml/Min) for water at 45 psig/3.1bar [membrane dia.]	
		13mm	25mm
0.2	42/2.8	22	175
0.45	30/2.1	NA	300

NA = Not available

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Luer-Lok is a registered trademark of Becton Dickinson. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.thermoscientific.com/filtration

ANZ: Australia: 1300 735 292, New Zealand: 0800 933 966;

Asia: China Toll-free: 800-810-5118 or 400-650-5118;

India: +91 22 6716 2200, India Toll-free: 1 800 22 8374;

Japan: +81-3-5826-1616; Other Asian countries: 65 68729717

Europe: Austria: +43 1 801 40 0; Belgium: +32 2 482 30 30;

Denmark: +45 4631 2000; France: +33 2 2803 2180;

Germany: +49 6184 90 6940, Germany Toll-free: 08001-536 376;

Italy: +39 02 95059 552 or 434-254-375;

Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 10200;

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 2000

8-0405-84 0212

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific