

# Thermo Scientific Nalgene Syringe Filters with 25mm Membranes

- Surfactant-Free Cellulose Acetate (SFCA)
- Cellulose Acetate (CA)
- Acrylic Housing

Cat. No.	Syringe Filter Description	Pore size, $\mu\text{m}$	Membrane Type	Packaging	No. per pack	No. per case
189-2000	Non-sterile prefilter	NA <sup>2</sup>	Glass-fiber	Bulk, zip-lock bag	—	50
190-2520 <sup>1</sup>	Sterile filter	0.2	SFCA	Individual blister pack	—	50
190-2545 <sup>1</sup>	Sterile filter	0.45	SFCA	Individual blister pack	—	50
190-2580 <sup>1</sup>	Sterile filter	0.8	CA	Individual blister pack	—	50
190-9920 <sup>1</sup>	Sterile filter	0.2	SFCA	Individual blister pack	—	125
190-9945 <sup>1</sup>	Sterile filter	0.45	SFCA	Individual blister pack	—	125
191-2020	Non-sterile filter	0.2	SFCA	Bulk, zip-lock bag	50	300
191-2045	Non-sterile filter	0.45	SFCA	Bulk, zip-lock bag	50	300
191-2080	Non-sterile filter	0.8	CA	Bulk, zip-lock bag	50	300
192-2520 <sup>1</sup>	Sterile prefilter/filter	NA <sup>2</sup> /0.2	CA	Individual blister pack	—	50

<sup>1</sup>Certified non-pyrogenic. <sup>2</sup>NA— not applicable.

**Thermo**  
SCIENTIFIC

Contact us for Sales and Service  
[thermoscientific.com/contactus](https://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

(For all syringe filters listed in this insert.)

1. Fill syringe with sample. (See "Caution")
2. Secure the syringe to the filter with a clockwise twist, then remove the syringe filter from its package.
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.

**NOTE:** To take advantage of the built-in glass-fiber prefilter on the upstream side of the CA membrane in Cat. No. 192, only pressure filtration should be used. DO NOT use aspiration or vacuum filtration with Cat. No. 192.

**Caution:** Use caution with syringes smaller than 10cc (10ml). The pressure generated in these syringes may exceed the 75-psig (5.1 bar) limit of the syringe filter.

### Aspiration or Vacuum Filtration

1. Secure the syringe to the filter with a clockwise twist, then remove the syringe filter from its package. 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.

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

Nalgene® syringe filters with cellulose acetate or surfactant-free cellulose acetate (CA or SFCA) membranes are specially designed to cold sterilize or filter small volumes of aqueous liquids using a syringe. CA and SFCA membranes offer low protein binding and are low in extractables. SFCA membranes have **no wetting agents or surfactants**. When used with a prefilter, overall filtration performance can be improved by up to 100%. They can be used with both standard slip-fit and Luer Lok® syringes and accept standard hypodermic needles.

Nalgene CA and SFCA membranes and syringe filters have good chemical resistance to aqueous solutions, dilute acids and bases.

They are **not recommended** for use with aldehydes, esters, alcohols, ketones or aliphatic, aromatic and halogenated

hydrocarbons or their derivatives. Refer to the current Nalgene Labware Catalog for further chemical resistance data.

Nalgene syringe filters are preassembled, disposable, non-cytotoxic and Triton™-free. Their unitary construction provides a leakproof seal and eliminates bypass. This minimizes fluid retention and will not compromise sterility. The acrylic filter housing supports a 25 mm membrane on both sides, allowing either pressure or vacuum filtration (aspiration). (*See note under Pressure Filtration.*)

**Syringe filters with CA or SFCA membranes are not autoclavable.**

Specifications					
Pore size (µm)	Max. oper. pressure (psig/bar)	Surface area, cm <sup>2</sup>	Bubble point (psig/bar)	Hold-up volume, ml	Nom. Flow rate for water at 14.7 psig/ 1 bar (ml/min)
0.2	75/5.1	5.3	45/3.1	<0.25	90
0.45	75/5.1	5.3	28/1.9	<0.25	180
0.8	75/5.1	5.3	7/0.5	<0.25	300

All syringe filters have a male Luer-Lok inlet and female Luer-Lok outlet.

©2009 Thermo Fisher Scientific Inc. All rights reserved. Luer Lok is a registered trademark of Becton-Dickinson. Triton is a trademark of Dow Chemical. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

[www.nalgenelabware.com](http://www.nalgenelabware.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 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 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

8-0403-81 0609

**Thermo**  
SCIENTIFIC