



## Thermo Scientific Nalgene Tank Liners

Ideal for single-use biopharmaceutical and diagnostic reagent fluid processing

- Designed specifically for Nalgene® Cylindrical Tanks
- Gamma-irradiated and non-pyrogenic
- USP Class VI, non-cytotoxic, food grade
- Sizes 19L - 757L (5 gal - 200 gal)

Thermo Scientific Nalgene Sterile Coex Polyethylene Film Tank Liners feature an open-bag, flat-bottom design to enhance mixing capabilities. The flexible liners offer cost-effective, single-use, biopharmaceutical and diagnostic reagent fluid processing using rigid Nalgene cylindrical plastic tanks. Coex Polyethylene film is free of animal-derived components (ADCF).

### Tank Liners, Sterile

Catalog No.	Liner Capacity	Fits Nalgene Tanks, Cat. No.	No. per case
343050-0005	19L / 5 Gal	11100-0005, 54100-0005	10
343050-0007	28L / 7.5 Gal	11100-0007, 54100-0007	10
343050-0010	38L / 10 Gal	11100-0010, 54100-0010	10
343050-0015	57L / 15 Gal	11100-0015, 54100-0015	10
343050-0030	113L / 30 Gal	11100-0030, 54100-0030	10
343050-0055	208L / 55 Gal	11100-0055, 54100-0055	10
343050-0080	303L / 80 Gal	11100-0080	10
343050-0100	378L / 100 Gal	11100-0100	10
343050-0150	568L / 150 Gal	11100-0150	10
343050-0200	757L / 200 Gal	11100-0200	10

10 individually heat-sealed packages in double polylined master carton for cleanliness.



© 2010 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

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

BRLSPNALGTANKLI 0610  
N20150 Ver. 1.0

**Thermo**  
SCIENTIFIC