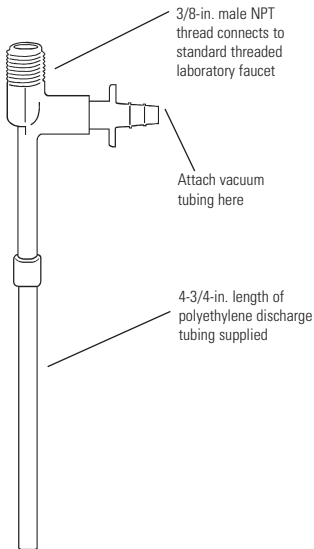


Thermo Scientific Nalgene Vacuum Pump (Aspirator type, Polypropylene)

Cat. No. 6140-0010

This Nalgene® vacuum pump has a maximum free air pumping capacity of 11.5 liters (3.0 gallons) per minute at a water flow rate of 6.5 liters (1.7 gallons) per minute. It will operate efficiently with water pressure as low as 7.5 psig (0.5 bar). The integral check valve eliminates vacuum loss after water is shut off. Maximum vacuum is 28.5 in. (724mm) Hg. To achieve maximum rate efficiency, use the tubing supplied with the pump and do not shorten the tubing length. If you cannot insert the tubing with moderate force and twisting action, hold the lower end of the pump under hot water for several seconds.



Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC

© 2010 Thermo Fisher Scientific Inc. All rights reserved. All 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 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

8-0400-29 0210

Thermo
SCIENTIFIC