

Thermo Scientific Nalgene Vacuum Plates

Cat.No. 5306-0070, 5306-0130

WARNING

Do not apply grease to either side of the neoprene gasket or to the lip of the Nalgene® vacuum jar.

The use of high vacuum grease, oil or any type of sealant ON EITHER SIDE of the gasket can result in movement and/or tearing of the gasket, the subsequent release of vacuum, ejection of the jar and possible injury.

After each use, check the gasket carefully for cuts or tears, and discard it if damaged.

For more information, contact Technical Support at:
technical.nalgene@thermofisher.com

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

www.thermoscientific.com

8-0402-26 1210

Thermo
SCIENTIFIC

Contact us for Sales and Service
thermoscientific.com/contactus

*Contact information contained within
this document may be incorrect.

Thermo
SCIENTIFIC