

# Thermo Scientific Nalgene Barbed Bulkhead Fittings

Cat. No. 6149-0001 (1/2 in. O.D.)

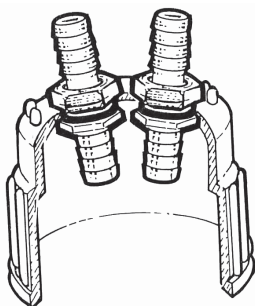
Cat. No. 6149-0002 (1/4 in. O.D.)

Nalgene® barbed bulkhead fittings permit the retrofitting of most Nalgene and other manufacturers' closures for liquid transfer when venting is not required.

Each kit includes:

- 2 barbed polypropylene fittings
- 2 acetal nuts
- 2 silicone gaskets
- 2 TPE (thermoplastic elastomer) port caps

All components are provided non-sterile. They may be sterilized by autoclave, gas or chemical methods.



## Installation Instructions

1. Determine the correct spacing of the holes by placing the applicable template on the selected closure.
2. For tubing with 1/2 in. I.D., use **Template A** (Cat. No. 6149-0001). Drill two 5/8 in. diameter holes approximately 1-1/2 inches apart in the top of the closure.
3. For tubing with 1/4 in. I.D., use **Template B** (Cat. No. 6149-0002). Drill two 3/8 in. diameter holes approximately 3/4 in. apart in the top of the closure.
4. Slide the gaskets past the threads on the fittings.
5. Push the fittings through the holes so that the gaskets are on the inside of the closure.
6. Screw the nuts onto the outside of the fittings.
7. Attach applicable tubing (not included). We recommend Nalgene 180 clear PVC plastic tubing or 50 silicone or 8060 silicone tubing.

## For vacuum applications

We recommend installing Nalgene barbed bulk head fittings on the Nalgene heavy-duty vacuum bottle (Cat. No. 2126-4000, -5000). This sturdy, heavy-walled, 4 or 5 liter bottle can withstand repeated applications of full vacuum for 24 hour periods at room temperature (20°C).

## For aseptic liquid transfer

Nalgene filling/venting closures, Cat. Nos. 2162-0830 and 2162-0831, fit all Nalgene carboys that accept an 83B closure. Supplied fully assembled, the filling/venting closures consist of a standard leakproof 83B Nalgene closure with TPE gasket and these modifications: two 1/2 in. or 1/4 in. barbed filling ports, one 3/8 in. vent port, two pieces of Nalgene 50 or 8060 silicone tubing and three TPE caps.

**Thermo**  
SCIENTIFIC

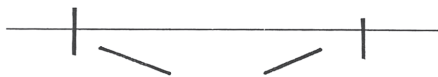
Contact us for Sales and Service  
[thermoscientific.com/contactus](http://thermoscientific.com/contactus)

\*Contact information contained within  
this document may be incorrect.

**Thermo**  
SCIENTIFIC

For further installation or application assistance, contact Technical Service at the numbers listed below.

## Template A



Drill 5/8 in. diameter holes here  
(centers are 1-1/2 in. apart)

## Template B



Drill 3/8-in. diameter holes here  
(centers are 3/4 in. apart)

© 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

8-0404-33 0310

**Thermo**  
SCIENTIFIC