



Thermo Scientific Nalgene Autoclaving Instructions

Recommended autoclave cycle is 121°C, 15 psig (1bar) for a minimum of 20 minutes. The specific cycle time needs to be determined and validated at each facility to assure sterility. Because plastics transfer heat more slowly than glass or metal, they will take longer to reach sterilization temperatures in the autoclave.

Carefully clean all items before autoclaving. This will prevent baking contaminants onto the surface of the plastic. We recommend Nalgene® L900 liquid detergent (Cat. No. 900-4000) for cleaning your plasticware. After cleaning, thoroughly rinse all items with distilled water.

Remove closures from all containers before autoclaving!

It is extremely important that the closure and container do not meet and seal, creating a vacuum inside the container. To prevent this (but keep the closure with its matching container), you should:

- Set the closure on top, but off-center of the container mouth.
- Use surgical wrap between the closure and container mouth to prevent threads from engaging.
- Tape the closure to the side of the container mouth.

Do not autoclave containers (except those made of fluorocarbons) containing detergent or wetting solutions. Do not autoclave products made of HDPE, LDPE, PETG, PS or PVC. Check for resin code on product.

After autoclaving, DO NOT screw on closure until the product has reached room temperature.











Printed on Rolland Enviro100 Print, which contains 100% post-consumer fiber, manufactured using renewable biogas energy and is certified EcoLogo, Processed Chlorine Free and FSC Recycled.

Visit www.thermoscientific.com product resources pages for complete product use guidelines.

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

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

www.thermoscientific.com

