



**NALGENE**

**Thermo**  
SCIENTIFIC

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

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

**This test tube rack is  
NOT AUTOCLAVABLE.**

[www.thermoscientific.com](https://www.thermoscientific.com)  
8-0405-85 0310

**Thermo**  
SCIENTIFIC